

Applied Biosystems SOLiD™ 4 System Templated Bead Preparation Guide

March 2010

Library
Preparation

**Templated Bead
Preparation**

Instrument
Operation



For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.

This user guide is the proprietary material of Applied Biosystems, LLC or its affiliates and is protected by laws of copyright. The customer of the SOLiD™ System is hereby granted limited, non-exclusive rights to use this user guide solely for the purpose of operating the SOLiD System. Unauthorized copying, renting, modifying, or creating derivatives of this user guide is prohibited.

Information in this document is subject to change without notice.

APPLIED BIOSYSTEMS DISCLAIMS ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE. TO THE FULLEST EXTENT ALLOWED BY LAW, IN NO EVENT SHALL APPLIED BIOSYSTEMS BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF, WHETHER OR NOT FORESEEABLE AND WHETHER OR NOT APPLIED BIOSYSTEMS IS ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

TRADEMARKS:

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

Covaris is a trademark of Covaris, Inc.

AmpliQ Gold and TaqMan are registered trademarks of Roche Molecular Systems, Inc. HydroShear is a registered trademark of Genomic Solutions, Inc. NanoDrop is a registered trademark of NanoDrop Technologies.

© Copyright 2010, Life Technologies Corporation. All rights reserved.

4448378 Rev. B

03/2010

Contents

	Preface	7
	Safety information	7
	How to use this guide	8
	How to obtain support	8
Chapter 1	Introduction	9
	Workflows	9
	Scales of preparation	12
Chapter 2	Templated Bead Preparation Overview	15
	Overview	16
	Organization of the protocol chapters	20
Chapter 3	Prepare Mini-Scale Templated Beads	21
	Prepare templated beads (mini-scale)	22
	Materials and equipment required (mini-scale)	22
	Workflow (mini-scale)	22
	Tips (mini-scale)	23
	Prepare the mini-scale emulsion PCR (ePCR) reaction	25
	Perform the emulsion break and bead wash (mini-scale)	33
	Enrich the mini-scale templated beads	37
	Modify the 3' ends (mini-scale)	41
Chapter 4	Prepare Full-Scale Templated Beads	45
	Prepare templated beads (full-scale)	46
	Materials and equipment required (full-scale)	46
	Workflow (full-scale)	46
	Tips (full-scale)	47
	Prepare the full-scale emulsion PCR (ePCR) reaction	49
	Perform the emulsion break and bead wash (full-scale)	57
	Enrich the full-scale templated beads	62
	Modify the 3' ends (full-scale)	66

Chapter 5	Macro-Scale Templated Beads (4 and 8 ePCR Reactions)	69
	Prepare templated beads (macro-scale)	70
	Materials and equipment required (macro-scale)	70
	Workflow (macro-scale)	71
	Tips (macro-scale)	72
	Macro-scale emulsion instructions	74
Chapter 6	Prepare Macro-Scale Templated Beads (4 ePCR Reactions)	75
	Prepare the templated beads (macro-scale: 4 ePCR reactions)	76
	Prepare 4 full-scale emulsion PCR (ePCR) reactions	76
	Perform the emulsion breaks and bead wash (macro-scale: 4 ePCR reactions)	76
	Enrich the templated beads (macro-scale: 4 ePCR reactions)	76
	Modify the 3' ends (macro-scale: 4 ePCR reactions)	82
Chapter 7	Prepare Macro-Scale Templated Beads (8 ePCR Reactions)	85
	Prepare the templated beads (macro-scale: 8 ePCR reactions)	86
	Prepare 4 full-scale emulsion PCR (ePCR) reactions	86
	Perform the emulsion breaks and bead wash (macro-scale: 8 ePCR reactions)	86
	Enrich the templated beads (macro-scale: 8 ePCR reactions)	86
	Modify the 3' ends (macro-scale: 8 ePCR reactions)	91
Appendix A	Required Materials	95
	Prepare templated beads (mini-scale)	96
	Prepare templated beads (full-scale and macro-scale)	100
Appendix B	Supplemental Procedures	105
	Program the Eppendorf Repeater® Xstream Pipettor	106
	Quantitate the beads using the NanoDrop® ND-1000 Spectrophotometer	107
	Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer	112
	Quantitate the beads using a hemocytometer	115
Appendix C	Library Concentration Conversion	117
Appendix D	Checklists and Workflow Tracking Forms	119
	Workflow checklists: prepare templated beads	120
	Workflow tracking: prepare templated beads (mini-scale or full-scale)	122
	Workflow tracking: prepare templated beads (macro-scale)	123

Appendix E	The Covaris™ S2 System	125
	Operation notes	126
	Covaris™ S2 programs	127
Appendix F	Instrument Warranty Information	129
	Computer configuration	130
	Limited product warranty	130
	Warranty period effective date	131
	Warranty claims	131
	Warranty exceptions	131
	Warranty limitations	132
	Damages, claims, and returns	133
Appendix G	Safety	135
	Instrumentation safety	136
	General instrument safety	136
	Physical hazard safety	136
	Chemical safety	137
	General chemical safety	137
	SDSs	138
	Chemical waste safety	139
	Biological hazard safety	141
	Glossary	143
	Documentation	147
	Related documentation	147
	Send us your comments	149
	Index	151

Safety information



Note: For important instrument safety information, refer to the *Applied Biosystems SOLiD™ 4 System Instrument Operation Guide* (PN 4448379). For general safety information, see this Preface and [Appendix G, “Safety” on page 135](#). When a hazard symbol and hazard type appear by a chemical name or instrument hazard, see [Appendix G, “Safety” on page 135](#) for the complete alert on the chemical or instrument.

Safety alert words

Four safety alert words appear in Applied Biosystems user documentation at points in the document where you need to be aware of relevant hazards. Each alert word—**IMPORTANT**, **CAUTION**, **WARNING**, **DANGER**—implies a particular level of observation or action, as defined below:



IMPORTANT! – Indicates information that is necessary for proper instrument operation, accurate chemistry kit use, or safe use of a chemical.



CAUTION! – Indicates a potentially hazardous situation that, if not avoided, may result in minor or moderate injury. It may also be used to alert against unsafe practices.



WARNING! – Indicates a potentially hazardous situation that, if not avoided, could result in death or serious injury.



DANGER! – Indicates an imminently hazardous situation that, if not avoided, results in death or serious injury. This signal word is to be limited to the most extreme situations.



SDSs

The Safety Data Sheets (SDSs) for any chemicals supplied by Applied Biosystems or Ambion are available to you free 24 hours a day. For instructions on obtaining SDSs, see [“SDSs” on page 138](#).



IMPORTANT! For the SDSs of chemicals not distributed by Applied Biosystems or Ambion contact the chemical manufacturer.

How to use this guide

- Text conventions** This guide uses the following conventions:
- **Bold** text indicates user action. For example:
Type **0**, then press **Enter** for each of the remaining fields.
 - *Italic* text indicates new or important words and is also used for emphasis.
For example:
Before analyzing, *always* prepare fresh matrix.
 - A right arrow symbol (▶) separates successive commands you select from a drop-down or shortcut menu. For example:
Select **File ▶ Open ▶ Spot Set**.
Right-click the sample row, then select **View Filter ▶ View All Runs**.
- User attention words** Two user attention words appear in Applied Biosystems user documentation. Each word implies a particular level of observation or action as described below:
-  **Note:** – Provides information that may be of interest or help but is not critical to the use of the product.
-  **IMPORTANT!** – Provides information that is necessary for proper instrument operation, accurate chemistry kit use, or safe use of a chemical.

How to obtain support

For the latest services and support information for all locations, go to:

www.appliedbiosystems.com

At the Applied Biosystems web site, you can:

- Access worldwide telephone and fax numbers to contact Applied Biosystems Technical Support and Sales facilities.
- Search through frequently asked questions (FAQs).
- Submit a question directly to Technical Support.
- Order Applied Biosystems user documents, SDSs, certificates of analysis, and other related documents.
- Download PDF documents.
- Obtain information about customer training.
- Download software updates and patches.

1

Introduction

Templated bead preparation is performed after library construction [refer to the *Applied Biosystems SOLiD™ 4 System Library Preparation Guide* (PN 4445673)]. To prepare templated beads, each library template is clonally amplified on SOLiD™ P1 DNA Beads by emulsion PCR (ePCR). After ePCR and enrichment of the templated beads, the templated beads are deposited onto a slide. The templates are then sequenced on the SOLiD™ 4 System [refer to the *Applied Biosystems SOLiD™ 4 System Instrument Operation Guide* (PN 4448379)].

Note: Customers who have purchased the SOLiD™ 4 EZ Bead™ System should not use this guide, but instead refer to the documentation delivered with the EZ Bead™ System.

Workflows

If you are preparing an ePCR reaction of a new library, you will obtain better sequencing results for a particular scale of templated bead preparation by titrating the library concentration to determine the optimal library concentration for ePCR (see [Figure 1 on page 10](#)).

To find the optimal library concentration, follow these steps:

1. Perform two separate ePCR reactions at library concentrations of 0.5 pM and 1.0 pM.
2. Perform a workflow analysis (WFA) run on the SOLiD™ 4 System to evaluate ePCR performance for each library concentration [refer to the *Applied Biosystems SOLiD™ 4 System Instrument Operation Guide* (PN 4448379)].
3. Use the optimal library concentration to prepare templated beads at the *same scale* of templated bead preparation you used to determine the optimal library concentration.

Note: You can determine the optimal library concentration for macro-scale templated bead preparation using the full-scale templated bead preparation.

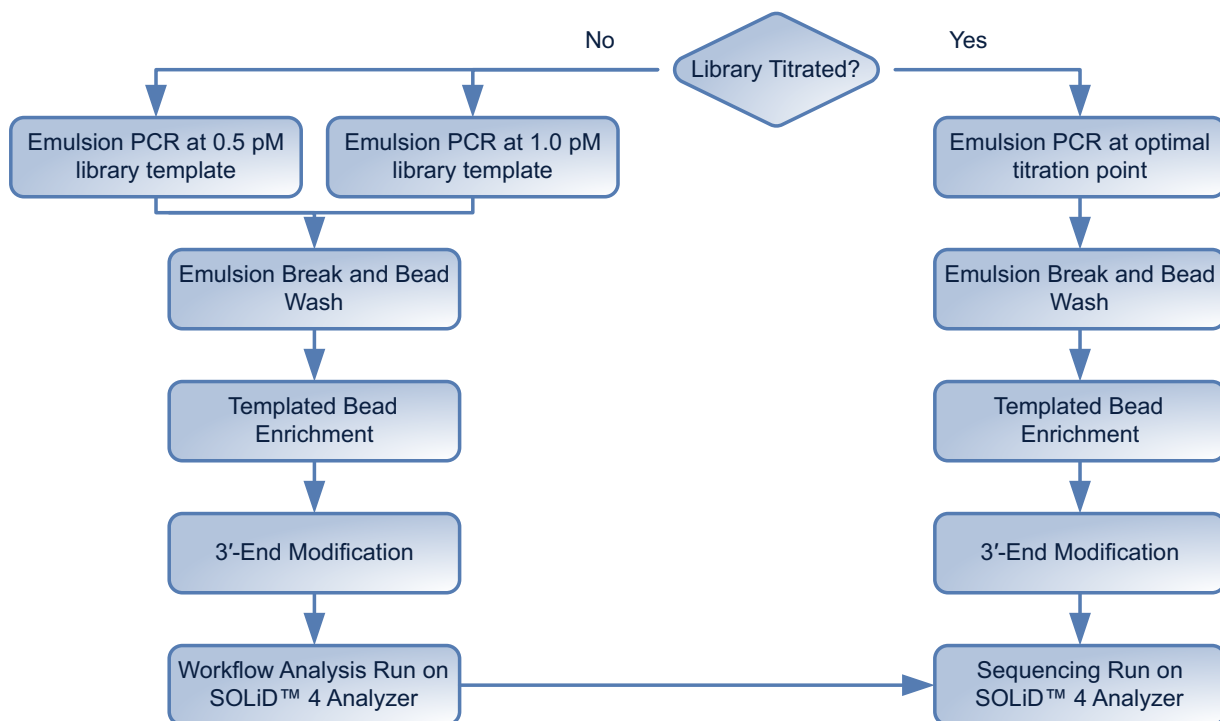


Figure 1 Workflow to prepare templated beads for SOLiD™ System sequencing using workflow analysis.

An alternative to preparing beads with two library titrations and performing a WFA run is quantitative PCR (qPCR). qPCR is a method to accurately measure library concentration. You can set up an emulsion PCR reaction according to the qPCR results because the molar optimal library concentration correlates with ePCR performance (see [Figure 2](#)). For details on qPCR, refer to the *Applied Biosystems SOLiD™ 4 System Library Preparation Guide* (PN 4445673). The workflow using qPCR is shown in the following figure.

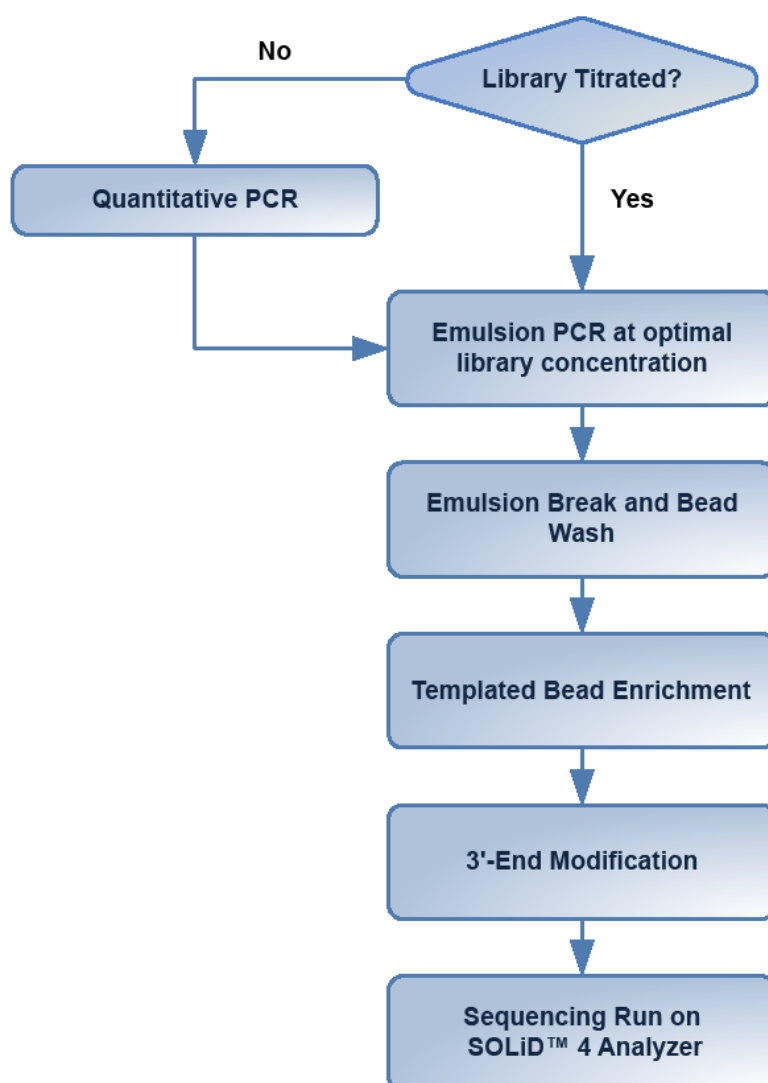


Figure 2 Workflow to prepare templated beads for SOLiD™ System sequencing using quantitative PCR.

Scales of preparation

You can prepare templated beads according to the amount of library that you want to amplify (see [Table 1](#)):

Table 1 Three ways to prepare templated beads according to the scale of preparation

Scale of preparation	Features	Go to ...
Mini	<ul style="list-style-type: none"> Yield: 75 to 150 million templated beads ePCR reaction: 1 ePCR reaction seeded with 800 million SOLiD™ P1 DNA Beads 	Chapter 3, Prepare Mini-Scale Templated Beads
Full	<ul style="list-style-type: none"> Yield: 150 to 300 million templated beads ePCR reaction: 1 ePCR reaction seeded with 1.6 billion SOLiD™ P1 DNA Beads 	Chapter 4, Prepare Full-Scale Templated Beads
Macro	<ul style="list-style-type: none"> Yield: 600 million to 1.2 billion templated beads ePCR reaction: 4 ePCR reactions, each seeded with 1.6 billion SOLiD™ P1 DNA Beads 	Chapter 5, Macro-Scale Templated Beads (4 and 8 ePCR Reactions) <i>and either</i> Chapter 6, Prepare Macro-Scale Templated Beads (4 ePCR Reactions)
	<ul style="list-style-type: none"> Yield: 1.2 billion to 2.4 billion templated beads ePCR reaction: 8 ePCR reactions, each seeded with 1.6 billion SOLiD™ P1 DNA Beads 	<i>or</i> Chapter 7, Prepare Macro-Scale Templated Beads (8 ePCR Reactions)

Choose the scale of templated bead preparation based on the number of templated beads required for the slide (see [Table 2](#)). Vary the targeted bead density for deposition based on your desired output, sample, and experimental conditions:

Table 2 Number of templated beads needed according to slide configuration

Slide configuration	Templated bead quantity requirements [‡] (millions per spot)
1-well	708
4-well	128
8-well	56

[‡] Assuming targeted bead deposition density of 300,000 beads per panel.

Decide which slide configuration is appropriate based on your desired output. Estimate expected output based on the number of beads, using the relationship shown below. Note that your actual output depends on your sample and the experimental conditions.

Expected output = (Number of templated beads) × (read length) × (% mappable beads)

Examples

1. For a fragment library with a 50-bp read length on 1 spot of an 8-well slide deposited at 300,000 beads/panel and assuming 50% matching:
Expected output = (56 million beads) \times (50 bp) \times 50% = 1.4 GB
2. For a mate-paired library with a 25-bp read length on a 1-well slide deposited at 300,000 beads/panel and assuming 60% matching:
Expected output = (708 million beads) \times (2 \times 25 bp) \times 60% = 21.2 GB

2

Templated Bead Preparation Overview

■ Overview..... 16

■ Organization of the protocol chapters..... 20

Overview

This guide describes how to clonally amplify short-fragment DNA populations onto SOLiD™ P1 DNA Beads using an emulsion method. Emulsions are made up of an oil phase containing emulsifiers and an aqueous phase, which includes PCR components (template, primers, DNA polymerase, and SOLiD™ P1 DNA Beads; see [Figure 3](#)).

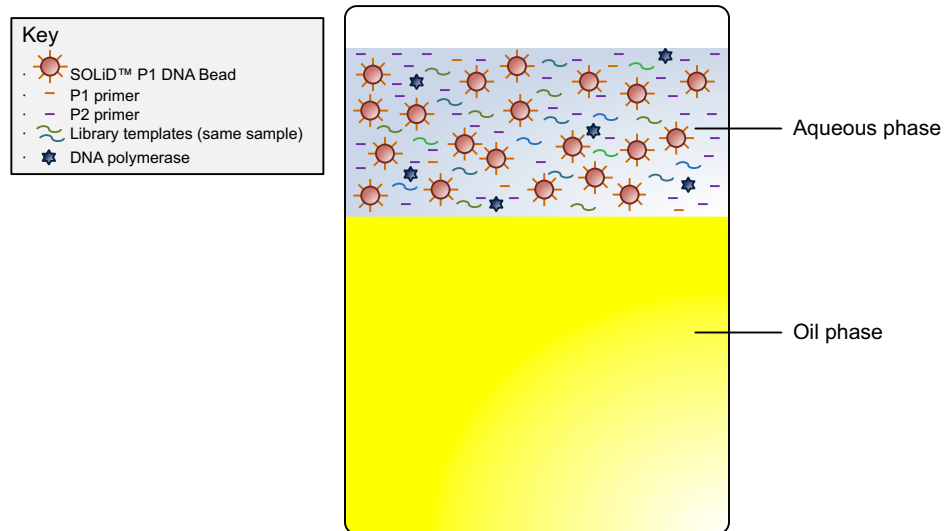


Figure 3 Aqueous phase and oil phase prior to the emulsification.

An emulsion is created using the ULTRA-TURRAX® Tube Drive from IKA®. An emulsion is made up of droplets of aqueous phase, or *micro-reactors*, in which the clonal amplification takes place. Micro-reactors containing a single SOLiD™ P1 DNA Bead and a single template, called *monoclonal micro-reactors*, are desired. However, Poisson bead distribution and Poisson template distribution allow for other types of reactors, including: *polyclonal micro-reactors* (contain that contain multiple templates); *non-clonal micro-reactors* (micro-reactors that contain no template); *multi-bead micro-reactors*; and micro-reactors with combinations of these characteristics (see [Figure 4](#) on page 17).

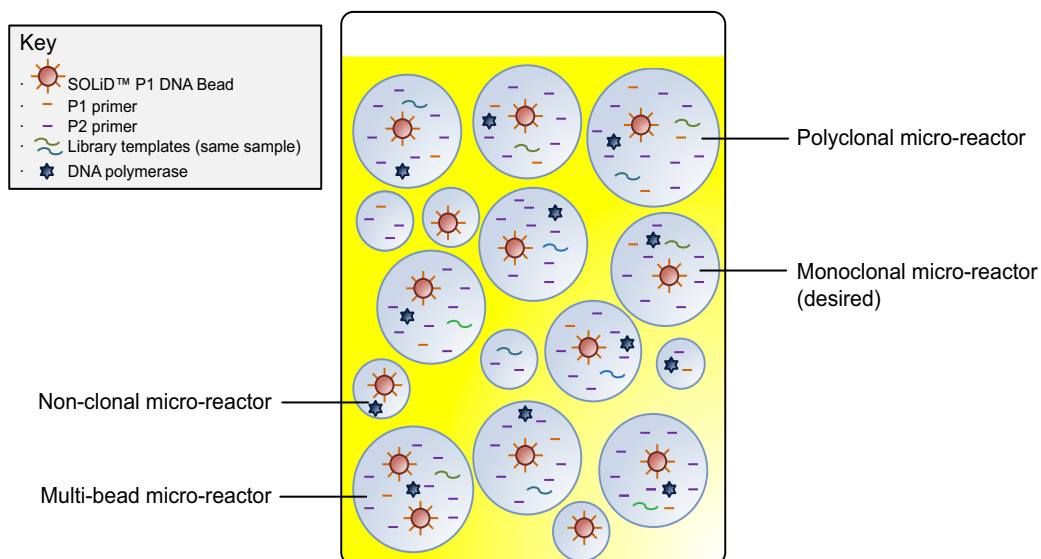


Figure 4 Emulsion before amplification (ePCR).

The emulsion is placed on a thermal cycler and run at standard PCR conditions. During emulsion PCR (ePCR), 30,000 or more copies of template are amplified onto each SOLiD™ P1 DNA Bead with the P1 Adaptor attached to the bead. In monoclonal and polyclonal micro-reactors, monoclonal and polyclonal templated beads are formed, respectively. In nonclonal micro-reactors, the SOLiD™ P1 DNA Bead cannot amplify. Multi-bead micro-reactors lead to suboptimal amplification (see [Figure 5](#)).

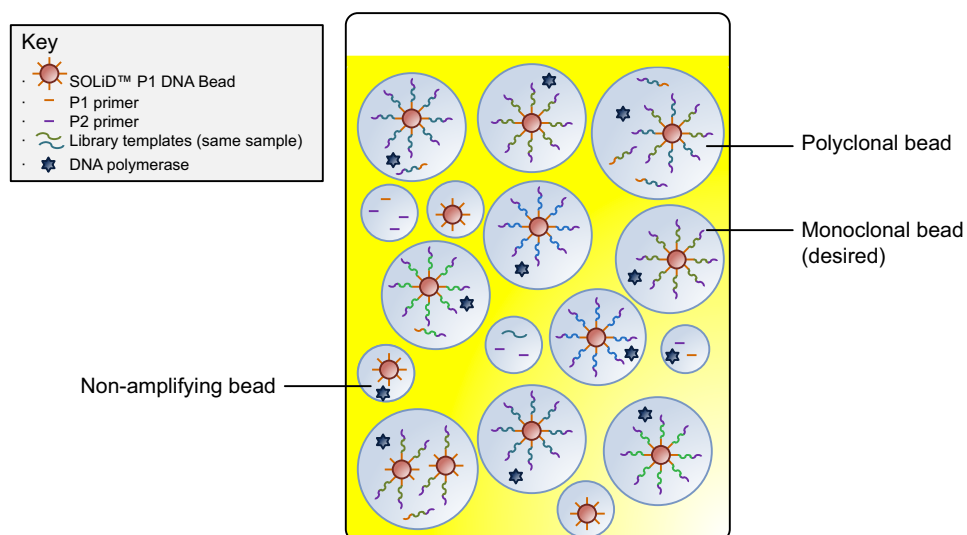


Figure 5 Emulsion after amplification (ePCR).

After emulsion PCR is complete, the micro-reactors in the emulsion are broken with 2-butanol, and the templated beads and nonamplifying beads are washed to clear away the oil and emulsifiers (see [Figure 6 on page 18](#)).

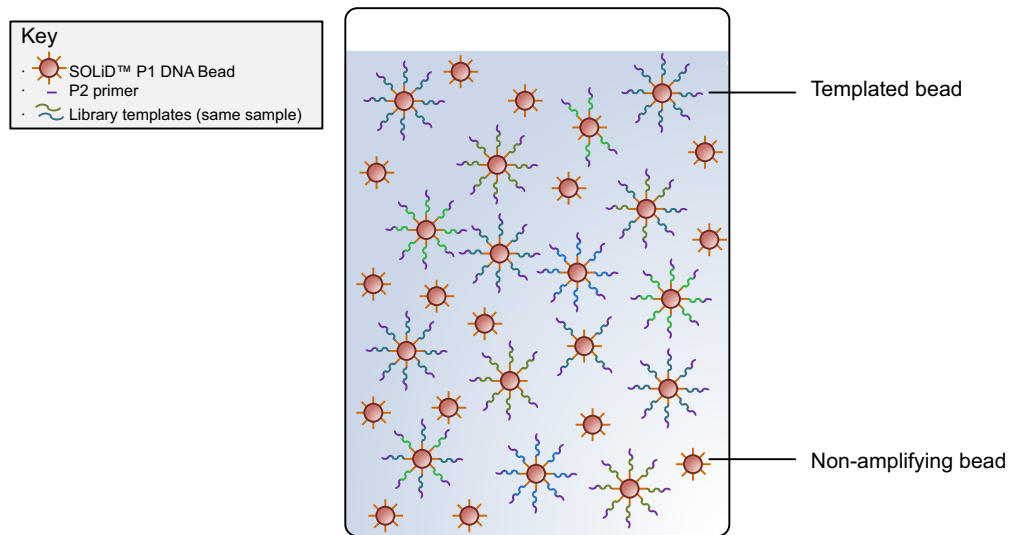


Figure 6 Templated and non-amplifying beads after emulsion break and bead wash.

Enrichment is required to isolate templated beads from non-amplifying or poorly amplifying beads. In an enrichment step, polystyrene beads with a single-stranded P2 Adaptor attached are used to capture templated beads. The mixture of enrichment beads, enrichment bead-templated bead complexes, and non-amplifying beads is centrifuged on a 60% glycerol cushion. The enrichment step results in a layer of enrichment beads (with or without templated beads attached) at the top and a layer of non-amplifying beads at the bottom. The layer of enrichment beads is extracted and denatured to dissociate the templated beads from the enrichment beads (see [Figure 7](#)).

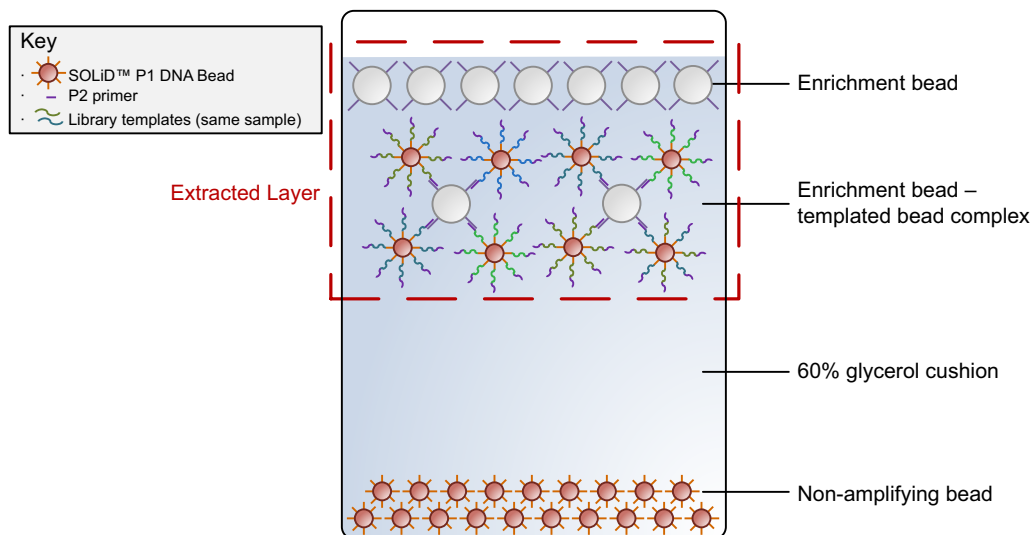


Figure 7 Enrichment beads and SOLiD™ P1 DNA beads after centrifugation with 60% glycerol.

In order to prepare the P2-enriched beads for deposition, a dUTP is added to the 3'-end of the P2 templates using a terminal transferase reaction (see [Figure 8](#)).

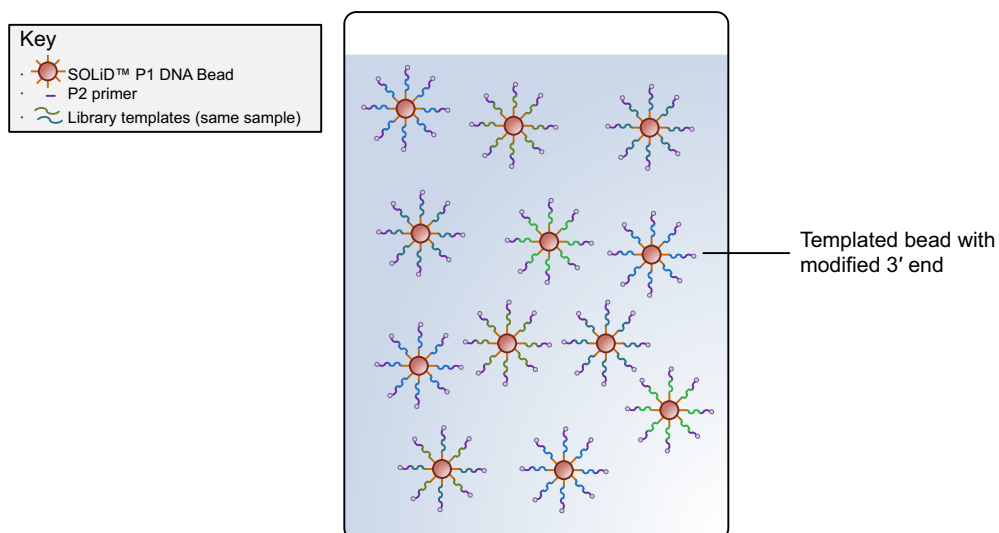


Figure 8 Templated beads after 3'-end modification.

Organization of the protocol chapters

The remaining chapters of this guide are organized by protocol scale:

Mini-Scale

[Chapter 3, “Prepare Mini-Scale Templated Beads” on page 21](#) describes how to generate 75 to 150 million templated beads using the *mini*-scale templated bead preparation method.

Full-Scale

[Chapter 4, “Prepare Full-Scale Templated Beads” on page 45](#) describes how to generate 150 to 300 million templated beads using the *full*-scale templated bead preparation method.

Macro-Scale (4 and 8 ePCR reactions)

[Chapter 5, “Macro-Scale Templated Beads \(4 and 8 ePCR Reactions\)” on page 69](#) provides a workflow overview and pre-emulsion PCR instructions for the *macro*-scale templated bead preparation method, which generates 600 million to 2.4 billion templated beads.

This chapter applies to macro-scale templated bead preparation with both 4 ePCR reactions and 8 ePCR reactions. To use the macro-scale templated bead preparation method, you must complete the instructions in [Chapter 5](#) before starting the instructions in either [Chapter 6](#) or [Chapter 7](#).

Macro-Scale (4 ePCR reactions)

[Chapter 6, “Prepare Macro-Scale Templated Beads \(4 ePCR Reactions\)” on page 75](#) describes how to generate 600 million to 2.4 billion templated beads using the *macro*-scale (4 ePCR reactions) templated bead preparation method.

The instructions in [Chapter 5](#) must be completed before you follow the instructions in [Chapter 6](#).

Macro-Scale (8 ePCR reactions)

[Chapter 7, “Prepare Macro-Scale Templated Beads \(8 ePCR Reactions\)” on page 85](#) describes how to generate 600 million to 2.4 billion templated beads using the *macro*-scale (8 ePCR reactions) templated bead preparation method.

The instructions in [Chapter 5](#) must be completed before you follow the instructions in [Chapter 7](#).

3

Prepare Mini-Scale Templated Beads

■ Prepare templated beads (mini-scale)	22
■ Materials and equipment required (mini-scale)	22
■ Workflow (mini-scale)	22
■ Tips (mini-scale)	23
■ Prepare the mini-scale emulsion PCR (ePCR) reaction	25
■ Perform the emulsion break and bead wash (mini-scale)	33
■ Enrich the mini-scale templated beads	37
■ Modify the 3' ends (mini-scale)	41

Prepare templated beads (mini-scale)

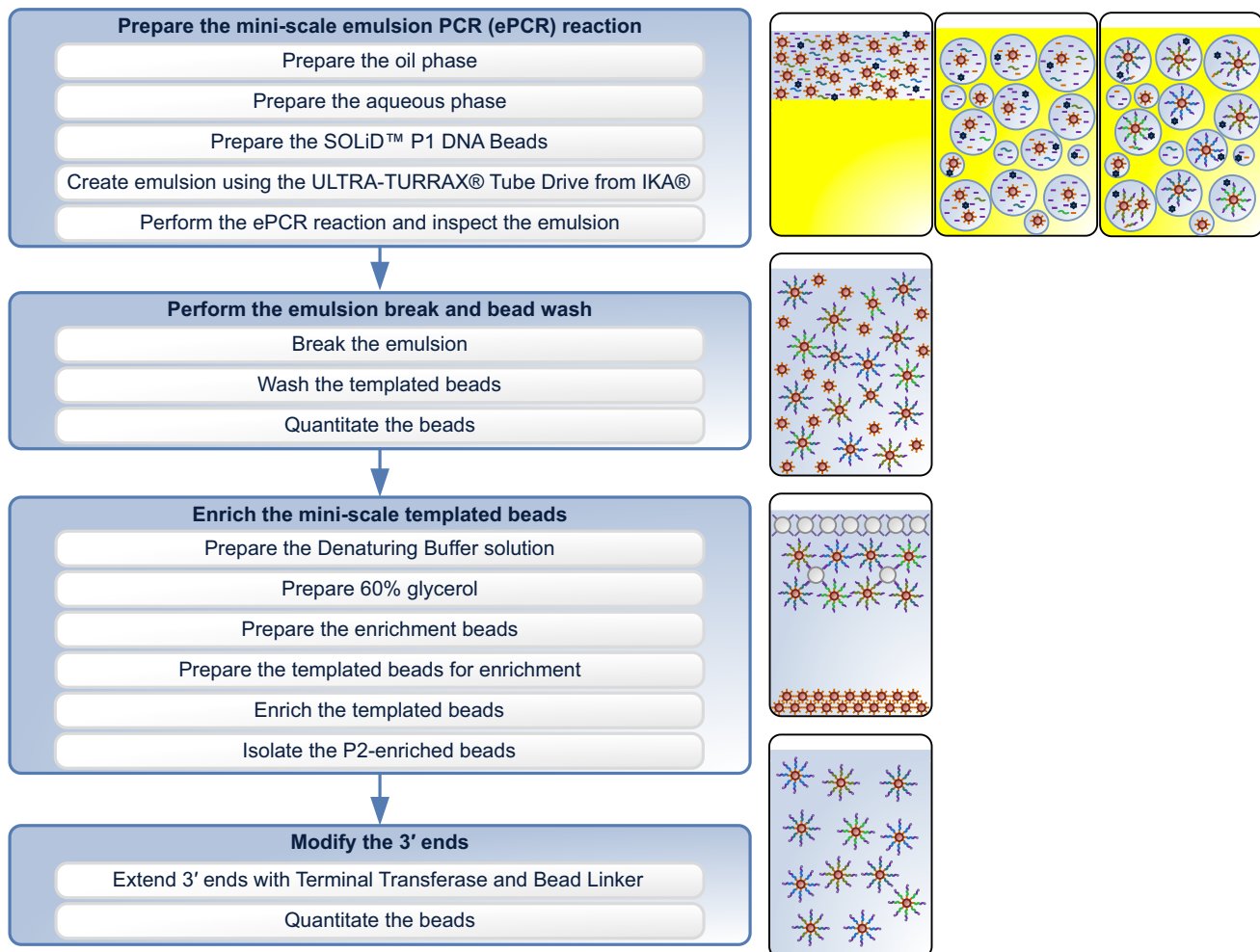
This chapter describes how to generate 75 to 150 million templated beads using the *mini-scale* templated bead preparation method.

Materials and equipment required (mini-scale)

See [Appendix A on page 95](#) for a list of equipment, kits, and consumables necessary for this procedure.

Workflow (mini-scale)

See the overview descriptions of the steps below the workflow.



**Workflow overview
(mini-scale)****Prepare the mini-scale emulsion PCR (ePCR) reaction**

The oil phase and aqueous phase of the emulsion are prepared separately, then emulsified using the ULTRA-TURRAX® Tube Drive from IKA®. Each emulsion is seeded with 800 million SOLiD™ P1 DNA Beads, then transferred into a single, 96-well plate for cycling. Different library template lengths require different numbers of cycles for thermal cycling.

Perform the emulsion break and bead wash (mini-scale)

The emulsion break uses 2-butanol to purify emulsified templated beads from the oil phase following amplification. The beads are washed to remove any residual 2-butanol, oil, and aqueous phase containing PCR reagents. The SOLiD™ Emulsion Collection Tray is placed over the 96-well plate, then the plate is centrifuged. Centrifuging the plate forces the emulsion from each well to a single reservoir. After centrifugation, 2-butanol is added to the reservoir. The broken emulsion is transferred to a 50-mL tube for further processing.

Enrich the mini-scale templated beads

The templated bead enrichment procedure isolates beads with full-length extension products following ePCR. Beads with full-length extension products are isolated by oligo hybridization using the sequence of the P2 primer. Both monoclonal and polyclonal beads are enriched. The procedure is designed to enrich for templated beads derived from one ePCR reaction yielding 75 to 150 million templated beads.

Modify the 3' ends (mini-scale)

The P2-enriched beads are extended with a Bead Linker by Terminal Transferase.

Tips (mini-scale)**General**

- Use syringes to accurately measure viscous reagents. Aspirate the volume very slowly from the reagent bottle to ensure that no air bubbles are trapped within the syringe. As is the best practice, draw some reagent into the syringe, dispense the entire reagent back to the reagent bottle, then draw the correct volume of reagent. Measure the volume to the point where the plunger contacts the side of the syringe.
- Perform all steps requiring 0.5-mL, 1.5-mL, and 2.0-mL tubes with Eppendorf LoBind Tubes. LoBind tubing from other vendors may have a chemical coating that has an adverse effect on bead deposition.
- Adjust microcentrifuge speeds and times according to the g-forces specified in the protocols. Applied Biosystems recommends the Eppendorf 5417R tabletop microcentrifuge.

**SOLiD™ P1 DNA
Beads**

- Do not freeze SOLiD™ P1 DNA Beads or templated beads. Store the SOLiD™ P1 DNA Beads at 4 °C in 1× TEX Buffer.

- If beads remain in the original tube after transfer, you can use a small additional volume of the appropriate buffer to recover the remaining beads. Do not exceed a total volume of 1.3 mL for a 1.5-mL LoBind Tube.

Covaris™ S2 System

- Applied Biosystems optimizes the procedures for the Covaris™ S2 System. The Covaris™ S2 System must be specially adapted to prepare beads for the Applied Biosystems SOLiD™ 4 System. Do not use the Covaris S1 sonicator or an unadapted Covaris™ S2 System for bead preparation. For more information, contact an Applied Biosystems SOLiD™ System applications specialist.
- To achieve optimal sonication by the Covaris™ S2 System, follow these guidelines:
 - Ensure that the Covaris™ S2 System is degassed.
 - Ensure that no bubbles are present in the Covaris™ S2 System.
 - Ensure that the instrument and tube are properly aligned for appropriate sonication of beads.
 - Use the appropriate adaptor with the Covaris™ S2 System, as shown in [Table 3](#). Place the tube collar at the indicator line of the adaptor.

Table 3 Tubes and adaptors for use with the Covaris™ S2 System

Sample volume	Size of LoBind Tube	Size of tube adaptor
≤ 200 µL	0.5-mL	0.65-mL
200 µL - 600 µL	1.5-mL	1.5-mL
600 µL - 1.2 mL	2.0-mL	1.5-mL

Prepare the mini-scale emulsion PCR (ePCR) reaction

Prepare the oil phase (mini-scale)

1. Use a 3-mL syringe to dispense 1.8 mL of Emulsion Stabilizer 1 into the 50-mL conical tube.
2. Use a 1-mL syringe to dispense 400 μ L of Emulsion Stabilizer 2 very slowly into the 50-mL tube.
3. Pour the Emulsion Oil (approximately 37.8 mL) into the tube that has the Emulsion Stabilizer 1 and Emulsion Stabilizer 2 so that the final volume is 40 mL.
4. Cap the 50-mL tube, then vortex the mixture until all Emulsion Stabilizer 1 and Emulsion Stabilizer 2 are incorporated into the Emulsion oil.
5. Allow the mixture to degas for a minimum of 20 minutes while you prepare the aqueous phase (see “[Prepare the aqueous phase \(mini-scale\)](#)”). To degas, place the mixture in a conical tube rack and slightly unscrew the conical tube cap.
6. Prime a 10-mL syringe by drawing in about 2 mL of oil phase then dispensing it back into the tube.
7. Take off the cap of a new SOLiD™ ePCR Tube. Use the primed 10-mL syringe to dispense 9 mL of oil phase into the SOLiD™ ePCR Tube, then cap the tube.

STOPPING POINT. The oil phase may be stored at 4 °C for up to 2 months. Before using the stored oil phase, thoroughly vortex and degas the solution for 20 minutes.

Prepare the aqueous phase (mini-scale)

1. Dilute ePCR Primer 1 to prepare a 10- μ M working stock solution. For each ePCR reaction, add 2 μ L of ePCR Primer 1 to 18 μ L of 1 \times Low TE Buffer. Mix well.
2. Using only 1 \times Low TE Buffer and LoBind Tubes, prepare a dilution of the library template to a final concentration of 500 pM. Use [Table 4 on page 26](#) to convert the mass/volume concentration to molar concentration for each library (for calculation details, see “[Library Concentration Conversion](#)” on page 117). Dilute only enough template for the desired number of emulsions. If needed, perform a serial dilution of the library to accurately obtain the desired library concentration. For example, perform a 5 \times dilution from 50 nM to 10 nM, then perform a 20 \times dilution from 10 nM to 500 pM.

Table 4 Concentration conversions by library type

Library type	Average library length (bp)	Molar concentration (pM)	Mass/volume concentration
Fragment Library	~215	500	71 pg/μL
2 x 50 bp Mate-Paired Library	~300	500	99 pg/μL
2 x 25 bp Mate-Paired Library	~155	500	51 pg/μL
Barcoded Fragment Library	~255	500	84 pg/μL
Whole Transcriptome Library	~230	500	76 pg/μL
Small RNA Library	~123	500	41 pg/μL
Barcoded SAGE Library	~130	500	43 pg/μL

! **IMPORTANT!** Do not freeze-thaw dilutions of the library more than 3 to 4 times. Stock solutions and dilutions of libraries should be stored at – 20 °C at a concentration of 5 ng/μL or greater.

- Choose the appropriate library concentration, then prepare the aqueous phase by combining the following reagents in a Nalgene wide-mouth jar according to the table below (see [Table 5](#)). For workflow analysis, prepare the aqueous phase for library concentrations of 0.5 pM and 1.0 pM.

Table 5 Prepare the aqueous phase

		Library concentration		
Component	Final concentration [‡]	0.5 pM	1.0 pM	X pM
		Volume per reaction (μL) [§]		
10X PCR Buffer	1X	280	280	280
dNTP Mix (100 mM mix comprised of 25 mM each of dATP, dTTP, dCTP, dGTP)	14 mM (3.5 mM of each dNTP)	392	392	392
Magnesium Chloride (1 M)	25 mM	70	70	70
ePCR Primer 1 (10 μM working stock solution)	40 nM	11.2	11.2	11.2
ePCR Primer 2 (500 μM)	3 μM	16.8	16.8	16.8
Template (500 pM)	0.5 pM or 1.0 pM	2.8	5.6	X × 5.6
Nuclease-free Water	N/A	1647.2	1644.4	1650 – (X × 5.6)
AmpliTaq Gold® DNA Polymerase, UP (5 U/μL)	0.54 U/μL	300	300	300
Total	N/A	2720	2720	2720

[‡] The final concentration is based on a total volume of 2800 μL, which includes 2720 μL of liquid components and 80 μL of beads.

[§] Volumes are for a single IKA®-based ePCR reaction to fill a 96-well plate.

Prepare the SOLiD™ P1 DNA Beads (mini-scale)

4. Keep the aqueous phase on ice until ready to use.

1. Thoroughly vortex one tube of SOLiD™ P1 DNA Beads. Invert the tube at least once during vortexing to ensure that any beads stuck to the cap are washed down, then pulse-spin the tube.
2. Place the tube of beads in the magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
3. Resuspend the beads in 200 µL of Bead Block Solution. Vortex the solution to ensure that all beads are suspended, then pulse-spin the tube.

⚠ **IMPORTANT!** Keep the Bead Block Solution at 4 °C until ready for use.

4. Sonicate the beads using the Bead Block Declump program on the Covaris™ S2 System (for program conditions, see [“Bead Block Declump” on page 127](#)), then pulse-spin the tube.
5. Place the tube in a magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
6. Resuspend the beads in 200 µL of 1× TEX Buffer and vortex to ensure that all beads are suspended, then pulse-spin the beads.

Create the emulsion with the ULTRA-TURRAX® Tube Drive from IKA® (mini-scale)

1. Place the SOLiD™ ePCR Tube containing 9 mL of oil phase onto the ULTRA-TURRAX® device, then twist the tube to lock it into position (see [Figure 9](#)).

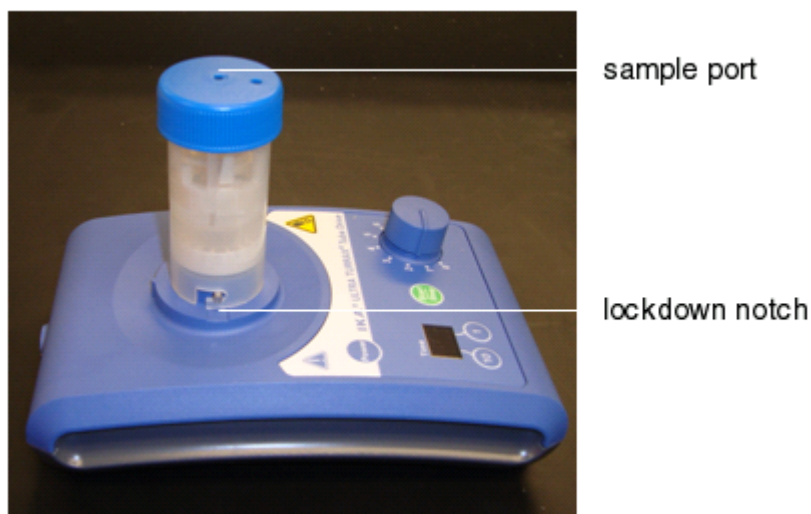


Figure 9 SOLiD™ ePCR Tube on the ULTRA-TURRAX® Tube Drive from IKA®.

2. Sonicate the SOLiD™ P1 DNA Beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)), then pulse-spin.

3. Immediately add 80 μL of the SOLiD™ P1 DNA Beads to the aqueous phase, then mix by gently swirling the bottle to ensure that the beads are uniformly dispersed (see [Figure 10](#)).



Figure 10 SOLiD™ P1 DNA Beads mixed in aqueous phase.

4. Verify that the Xstream pipettor is set up for mini-scale emulsions (see [Figure 11](#)):
 - Dial Setting: **Pip**
 - Speed (aspirate UP): **scale 5** (mid-range)
 - Speed (dispense DOWN): **scale 1** (slowest)
 - Total volume: **2.80 mL**

If necessary, reprogram the Xstream pipettor (see “[Program the Eppendorf Repeater® Xstream Pipettor](#)” on page 106).



Figure 11 Xstream pipettor settings, with the dial setting set to Pip.

5. Attach a 10-mL Combitip Plus tip onto the Xstream pipettor.
6. Fill the 10-mL Combitip Plus tip with the entire 2.80 mL of aqueous phase and bead mixture with the Xstream pipettor (see [Figure 12 on page 29](#)).



Figure 12 Filling the 10-mL Combipip Plus tip with the aqueous phase and bead mixture using the Xstream pipettor.

7. Verify that the time on the ULTRA-TURRAX[®] Tube Drive from IKA[®] is set to **5** minutes, then press the **Start** button.
8. Wait for the instrument's fly wheel to engage and to reach proper speed, then gently place the Combipip Plus tip into the center sample loading hole in the ULTRA-TURRAX[®] cap (see [Figure 13 on page 30](#)).



Figure 13 Correct placement of Combitip Plus into sample port in SOLiD™ ePCR Tube cap.

9. Dispense the aqueous phase and bead mixture into the spinning oil phase. When the entire volume is dispensed, press the center blue button *twice* on the pipettor to empty all contents from the Combitip Plus tip.
10. Remove a 5-mL Combitip Plus tip from its packaging, then cut off its end at the bevel with a razor blade (see [Figure 14](#)).

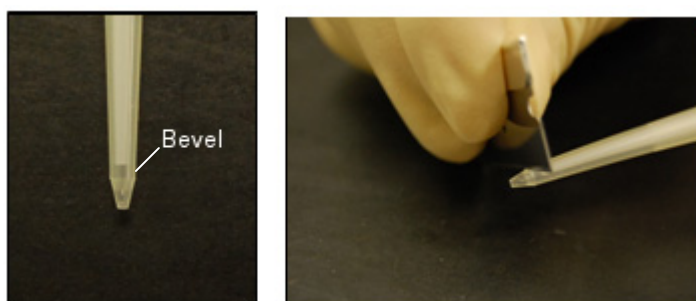


Figure 14 Cutting the Combitip Plus tip for emulsion dispersion.

11. Attach the cut Combitip Plus tip onto an Eppendorf Repeater® Plus Pipette.

12. Gently dispense 100 μ L of emulsion into each well of a 96-well PCR plate, then seal the plate with clear adhesive film (see [Figure 15](#)).



Figure 15 Emulsion transferred to a 96-well plate.

Perform the ePCR
reaction and
inspect the
emulsion (mini-
scale)

1. Set up the ePCR conditions on the GeneAmp[®] PCR System 9700:
 - ePCR thermal cycling program:

Stage	Step	Temp (°C)	Time
Holding	Denature	95	5 min
40 cycles [‡] or 60 cycles [§]	Denature	93	15 sec
	Anneal	62	30 sec
	Extend	72	75 sec
Holding	Final extension	72	7 min
Holding	—	4	∞

[‡] Set 40 cycles: Fragment library or 2 \times 25 bp mate-paired library.

[§] Set 60 cycles: 2 \times 50 bp mate-paired library.

- Ramp speed: **9600**
 - Reaction volume: **50 μ L**
2. Place the 96-well plate in a GeneAmp[®] PCR System 9700, then start the run.
 3. After the ePCR Program finishes, inspect the bottom of the reaction plate for beads that have fallen out of the emulsion. Beads appear as amber-colored specks at the bottom of a well. A small number of beads may fall out of emulsion and appear as small brown flecks at the bottom of a well. Applied Biosystems does not recommend any further processing of emulsions that have more than 3 wells of broken emulsion, where aqueous phase appears at the bottom of a well (see [Figure 16 on page 32](#)).

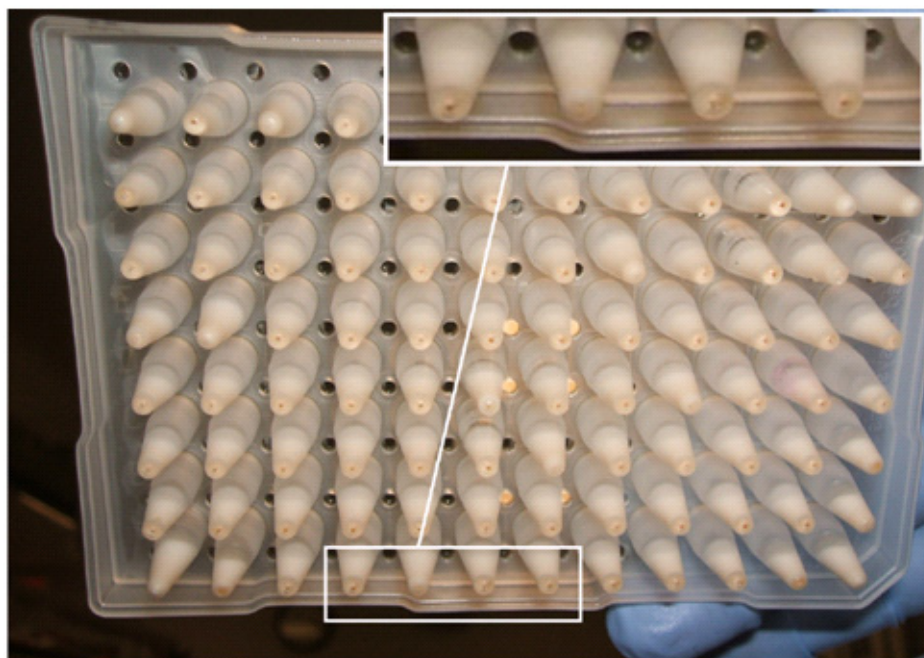


Figure 16 Broken emulsions.

STOPPING POINT. Store the 96-well plate at 4 °C, or proceed to [“Perform the emulsion break and bead wash \(mini-scale\)”](#) on page 33.

Perform the emulsion break and bead wash (mini-scale)

Break the emulsion (mini-scale)

1. Place the SOLiD™ Emulsion Collection Tray on top of the ePCR 96-well plate (see [Figure 17](#)).

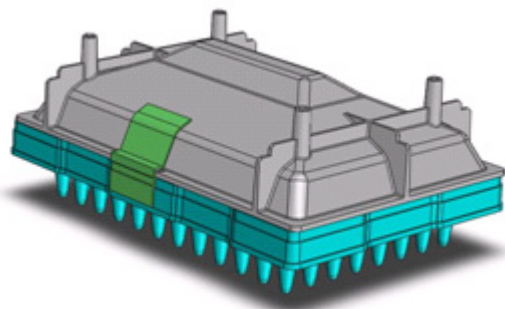


Figure 17 The SOLiD™ Emulsion Collection Tray taped to a 96-well reaction plate.

2. Seal the pieces together with tape on all four sides. Immediately prior to centrifugation, flip the entire apparatus so that the 96-well plate is upside-down over the collection tray (see [Figure 18](#)).

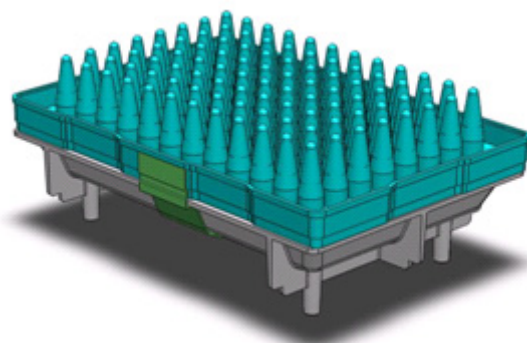


Figure 18 The inverted plate assembly.

3. Centrifuge the inverted plate and reservoir for 2 minutes at $550 \times g$ according to the recommended centrifuge settings (see [Table 6](#)).

Table 6 Recommended centrifuge settings

Adjustable Parameter	Recommend Setting
Acceleration (independent)	High
Deceleration (independent)	Low
Acceleration/Deceleration (single setting)	Mid

4. After centrifugation is complete, remove the plate assembly from the centrifuge and place on the lab bench. Hold the assembly steady, then gently remove the tape and 96-well plate from the collection tray.



Note: Ensure that the centrifuge is working properly and maintained regularly. Use anti-slip pads in the centrifuge carriers whenever possible.

5. In a fume hood, add 10 mL of 2-butanol to the collection tray using a serological pipette.
6. Pipet the emulsion up and down until the solution is homogeneous.
7. Transfer all the emulsion and 2-butanol to a 50-mL conical tube.
8. Rinse the reservoir with an additional 6 mL of 2-butanol to ensure that all residual beads are recovered.
9. Cap the tube, then vortex to mix the solution.
10. Centrifuge the tube at $2000 \times g$ for 5 minutes.
11. Gently decant the 2-butanol-oil phase into a waste bottle. With the tube inverted, place the tube onto paper towels to drain residual 2-butanol oil.
12. Wait 5 minutes to ensure that all the oil is removed.

Wash the templated beads (mini-scale)

1. Place the 50-mL tube upright in a rack, then add 600 μ L of 1 \times Bead Wash Buffer. Let the pellet soak in 1 \times Bead Wash Buffer for 2 minutes.
2. Resuspend the pellet by gently pipetting up and down, then transfer the beads from the 50-mL tube to a 1.5-mL LoBind Tube.
3. Rinse the *bottom* of the 50-mL tube with an additional 600 μ L of 1 \times Bead Wash Buffer, then transfer the wash to the 1.5-mL LoBind Tube.
4. Vortex the 1.5-mL LoBind Tube, then centrifuge the tube at $21,000 \times g$ (minimum $14,000 \times g$) for 1 minute.
5. Remove the top oil phase with a pipette. Remove as much of the oil at the meniscus as possible.
6. With a new pipette tip, carefully remove and discard the supernatant.
7. Resuspend the pellet by adding 150 μ L of 1 \times Bead Wash Buffer to the tube, then vortex the tube. Pulse-spin the tube, then transfer the mixture to a new 1.5-mL LoBind Tube.
8. Rinse the *bottom* of the original tube with an additional 150 μ L of 1 \times Bead Wash Buffer, then transfer the wash to the new tube.

9. Add 1 mL of 1× Bead Wash Buffer to the new tube, the vortex the tube.
10. Centrifuge the tube at $21,000 \times g$ (minimum $14,000 \times g$) for 1 minute, then remove and discard the supernatant.
11. Resuspend the beads in 200 μL of 1× TEX Buffer.
12. Place the tube in a magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
13. Resuspend the beads in 200 μL of 1× TEX Buffer.

STOPPING POINT. Store the templated beads at 4 °C in 1× TEX Buffer, or proceed to “Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer (mini-scale)” or “Quantitate the beads using a hemocytometer” on page 115.

Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer (mini-scale)

1. If necessary, generate a standard curve (see “Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer” on page 112).
2. Sonicate the beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see “Covalent Declump 1” on page 127), then pulse-spin the beads.
3. Make a 1-mL dilution of beads in 1× TEX Buffer (1:10 dilution recommended) in a 1.5-mL LoBind Tube.
4. Use the SOLiD™ Bead Concentration Chart (Applied Biosystems PN 4415131) to estimate the bead concentration of the beads (see Figure 19 for a picture of the chart; for best results use the official chart).

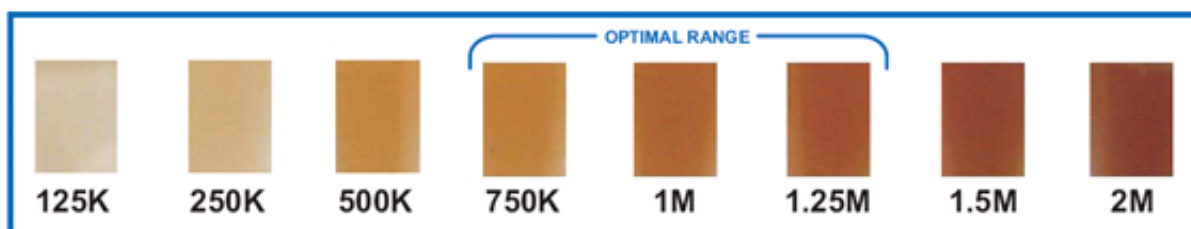


Figure 19 The SOLiD™ Bead Concentration Chart (facsimile). For best results, use the SOLiD™ Bead Concentration Chart (PN 4415131), supplied separately.

5. Adjust the volume of beads so that the color of the bead solution matches a color in the optimal range (750,000 beads/ μL to 1.25 million beads/ μL). See Figure 20 on page 36 for the workflow.

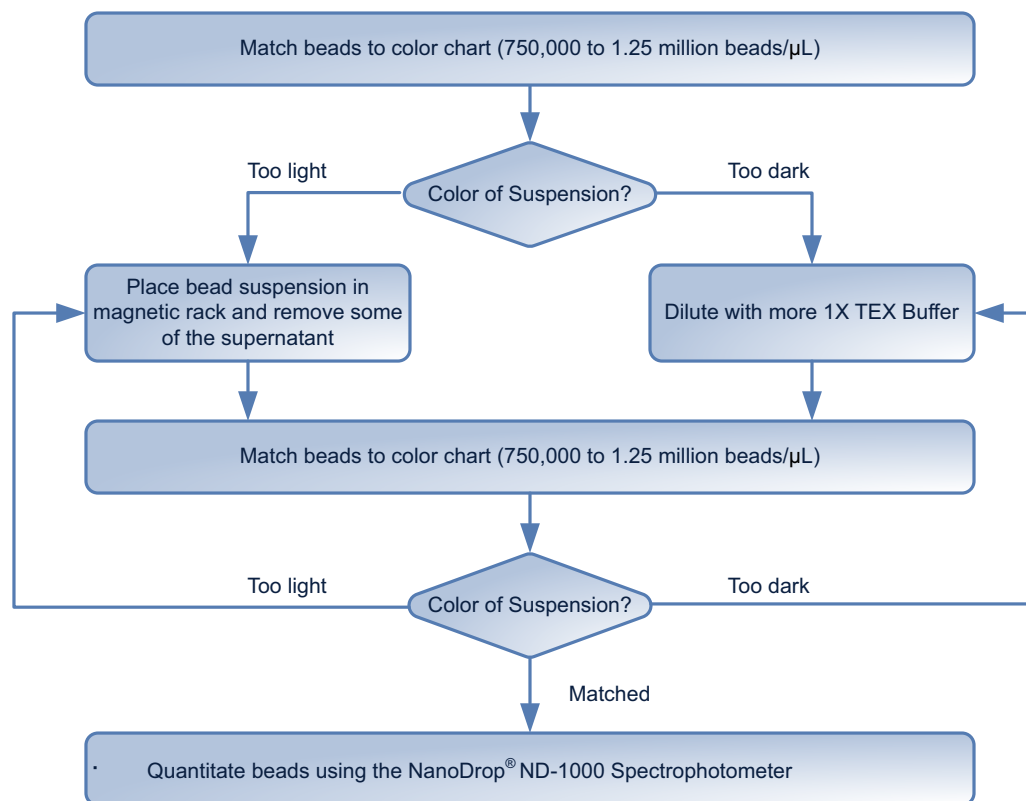


Figure 20 The SOLiD™ Bead Concentration Chart workflow.

6. When the bead concentration is within accurate range, quantitate the beads using the NanoDrop® ND-1000 Spectrophotometer. Take three readings, then average them. Calculate the bead concentration using the appropriate standard curve (for more details, see [“Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer”](#) on page 112).
7. Combine all diluted and undiluted beads.

STOPPING POINT. Store the beads at 4 °C in 1X TEX Buffer, or proceed to [“Enrich the mini-scale templated beads”](#) on page 37.

Enrich the mini-scale templated beads

Prepare the Denaturing Buffer solution (mini-scale)

1. For each ePCR reaction, transfer 1.8 mL of Denaturing Buffer to a 15-mL conical tube.
2. Add 200 μ L of Denaturant to the 1.8 mL of Denaturing Buffer, then cap the tube and vortex.

⚠ **IMPORTANT!** Prepare a new batch of the prepared Denaturing Buffer solution each week.

Prepare 60% glycerol (mini-scale)

1. With a 10-mL syringe, add 4 mL of Nuclease-free Water to a 15-mL conical tube.
2. With a 3-mL syringe, add 6 mL of glycerol to the Nuclease-free Water by dispensing 3 mL of glycerol twice with the syringe. Fill and dispense the glycerol slowly to ensure that the total volume of glycerol is dispensed.
3. Cap the tube, then vortex to mix the solution well.

⚠ **IMPORTANT!** Prepare the 60% glycerol fresh weekly.

Prepare the enrichment beads (mini-scale)

1. Vortex the enrichment beads and immediately transfer 300 μ L of the enrichment beads to a 1.5-mL LoBind Tube.
2. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
3. Resuspend the enrichment beads in 900 μ L of 1 \times Bind & Wash Buffer.
4. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
5. Repeat steps 3 and 4.
6. Resuspend the enrichment beads in 150 μ L of 1 \times Bind & Wash Buffer.
7. Add 1.5 μ L of 1 mM Enrichment Oligo to the tube of enrichment beads, then vortex and pulse-spin the tube.
8. Rotate the tube at room temperature for 30 minutes.
9. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
10. Resuspend the beads in 900 μ L of 1 \times TEX Buffer.

11. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
12. Repeat steps 10 and 11.
13. Resuspend the enrichment beads in 75 μL of 1 \times Low Salt Binding Buffer.

STOPPING POINT. Store the prepared enrichment beads at 4 °C in 1 \times Low Salt Binding Buffer, or proceed to [“Prepare the templated beads for enrichment \(mini-scale\)”](#). Prepared enrichment beads should be used within one week of preparation

Prepare the templated beads for enrichment (mini-scale)

1. Place the tube in the magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
2. Resuspend the templated beads in 300 μL of prepared Denaturing Buffer solution, then let the suspension stand for 1 minute.
3. Place the tube in the magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
4. Repeat steps 2 and 3 twice.
5. Resuspend the templated beads in 300 μL of 1 \times TEX Buffer.
6. Place the tube in the magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
7. Repeat steps 5 and 6 twice.
8. Resuspend the templated beads in 75 μL of 1 \times TEX Buffer, then transfer the templated bead suspension to a new 0.5-mL LoBind Tube.
9. Sonicate the beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)).

Enrich the templated beads (mini-scale)

1. Transfer the prepared enrichment beads to the tube of templated beads, then vortex and pulse-spin the bead mixture.
2. Sonicate the enrichment-templated bead mixture using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 3” on page 127](#)), then pulse-spin the beads.
3. Incubate the bead mixture at 61 °C for 15 minutes. During the incubation, vortex and pulse-spin the bead mixture every 5 minutes including at the end of the incubation.
4. Immediately cool the beads on ice for 2 minutes.
5. Add 400 μL of *freshly prepared* 60% glycerol to a new 1.5-mL LoBind Tube.

6. Gently pipet the bead mixture up and down to mix, then load the entire volume *carefully* on top of the 60% glycerol solution. Do *not* vortex the tube.
7. Centrifuge the tube for 3 minutes at $21,000 \times g$ (minimum $14,000 \times g$).
8. Add 1 mL of 1× TEX Buffer to a new 2.0-mL LoBind Tube.
9. Transfer the top layer of beads into the tube with 1× TEX Buffer. Aspirate as little glycerol as possible to collect all of the beads at the top layer without touching the un-templated beads at the bottom of the tube. When you dispense the top layer of beads into the 1× TEX Buffer, dispense the beads into the bottom of the tube. Aspirate a small amount of 1× TEX buffer to clean the pipette tip.
10. Top off the tube with additional 1× TEX Buffer to the 2.0-mL mark, then vortex.
11. Centrifuge the tube for 1 minute at $21,000 \times g$ (minimum $14,000 \times g$).



Note: Verify that the beads are pelleted. Excess glycerol carried over to the 1× TEX Buffer creates a matrix that impedes pelleting of beads.

Proceed according to the table below (see [Table 7](#)):

Table 7 Steps for pelleted or unpelleted beads

If the beads are...	Then perform steps...
Pelleted	12 and 13
Not pelleted	14 to 16

12. Remove the supernatant. Add 400 μ L of 1× TEX Buffer to the tube of beads and vortex.
13. Proceed to [“Isolate the P2-enriched beads \(mini-scale\)” on page 40](#).
14. Transfer half of the tube volume to a new 2.0-mL LoBind Tube, then add an additional 500 μ L of 1× TEX Buffer to each tube. Vortex each tube.
15. Centrifuge the tubes for 1 minute at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
16. Add 200 μ L of 1× TEX Buffer to each tube, resuspend the beads, then pool the beads into one tube.

**Isolate the
P2-enriched beads
(mini-scale)**

1. Centrifuge the tube for 1 minute at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.

❗ **IMPORTANT!** *Never* magnet the P2-enriched beads before adding prepared Denaturing Buffer solution to the beads. If you do, the templated beads linked to the enrichment beads are lost when the supernatant is removed.
2. Resuspend the pellet with 400 μL of prepared Denaturing Buffer solution, then let the solution stand for 1 minute.
3. Place the tube in a magnetic rack for at least 1 minute until the supernatant is pure white or clear, then remove and discard the supernatant.
4. Repeat steps 2 and 3 until the supernatant is clear (all white enrichment beads have been removed).
5. Resuspend the beads in 400 μL of 1 \times TEX Buffer.
6. Place the tube in a magnetic rack for at least 1 minute until the supernatant is clear. Remove and discard the supernatant.
7. Repeat steps 5 and 6.
8. Resuspend the beads in 200 μL of 1 \times TEX Buffer. Vortex, pulse-spin, then transfer the bead solution to a 1.5-mL LoBind Tube.
9. Rinse the 2.0-mL tube with 200 μL of 1 \times TEX Buffer and transfer the rinse to the 1.5-mL LoBind Tube.
10. Sonicate the beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)). Pulse-spin the beads.
11. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
12. Resuspend the beads in 400 μL of 1 \times TEX Buffer.
13. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
14. If the supernatant appears cloudy due to residual enrichment beads, repeat steps 12 and 13 until the supernatant is clear.
15. Resuspend the beads in 400 μL of 1 \times TEX Buffer.

STOPPING POINT. Store the prepared enrichment beads at 4 °C in 1 \times TEX Buffer, or proceed to [“Modify the 3’ ends \(mini-scale\)”](#).

Modify the 3' ends (mini-scale)

Extend the 3' ends with Terminal Transferase and Bead Linker (mini- scale)

1. If the P2-enriched beads have been stored overnight or longer, sonicate the beads using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 3” on page 127](#)). Pulse-spin the beads.
2. For each ePCR reaction, prepare 500 µL of 1× Terminal Transferase Reaction Buffer (see [Table 8](#)):

Table 8 Prepare Terminal Transferase Reaction Buffer

Component	Volume per reaction (µL)
10× Terminal Transferase Buffer	55
10× Cobalt Chloride	55
Nuclease-free Water	390
Total	500



Note: The 1× Terminal Transferase Reaction Buffer should be clear. If the solution becomes colored, discard it and prepare a fresh buffer using a new lot of material.

3. Add 1 µL of 50 mM Bead Linker to 49 µL of 1× Low TE Buffer to prepare a 1 mM Bead Linker solution.
4. Place the tube of P2-enriched beads in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
5. Resuspend the beads in 100 µL of 1× Terminal Transferase Reaction Buffer, then transfer the beads to a new 1.5-mL LoBind Tube.
6. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
7. Resuspend the beads in 100 µL of 1× Terminal Transferase Reaction Buffer.
8. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
9. Resuspend the beads in 178 µL of 1× Terminal Transferase Reaction Buffer.
10. Add 20 µL of 1 mM Bead Linker solution.
11. Sonicate the beads using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 3” on page 127](#)). Pulse-spin the beads.
12. Add 2 µL of Terminal Transferase (20 U/µL) and vortex. Pulse-spin the beads.

13. Place the tube on a rotator and rotate for 2 hours at 37 °C.
14. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
15. Resuspend the beads in 400 µL of 1× TEX Buffer.
16. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
17. Resuspend the beads in 200 µL of 1× TEX Buffer.

STOPPING POINT. Store the templated beads at 4 °C in 1× TEX Buffer, or proceed to [“Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer \(mini-scale\)”](#) or [“Quantitate the beads using a hemocytometer”](#) on page 115.

Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer (mini-scale)

1. If necessary, generate a standard curve (see [“Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer”](#) on page 112).
2. Sonicate the beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1”](#) on page 127), then pulse-spin the beads.
3. Use the SOLiD™ Bead Concentration Chart (Applied Biosystems PN 4415131) to estimate the bead concentration of the beads (see [Figure 21](#) for a picture of the chart; for best results use the official chart).

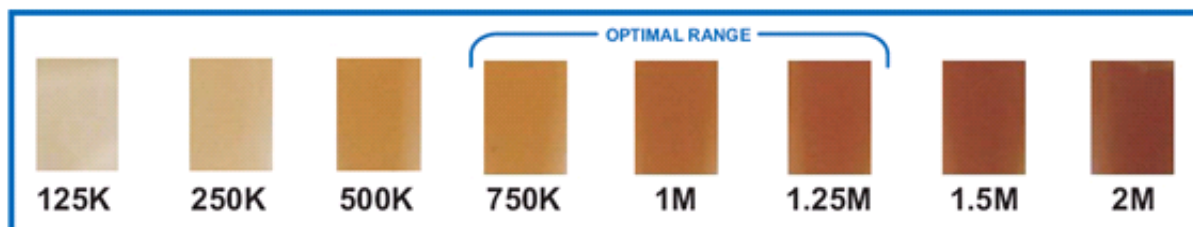


Figure 21 The SOLiD™ Bead Concentration Chart (facsimile). For best results, use the SOLiD™ Bead Concentration Chart (PN 4415131), supplied separately.

4. Adjust the volume of beads so that the color of the bead solution matches a color in the optimal range (750,000 beads/µL to 1.25 million beads/µL). See [Figure 22 on page 43](#) for the workflow.

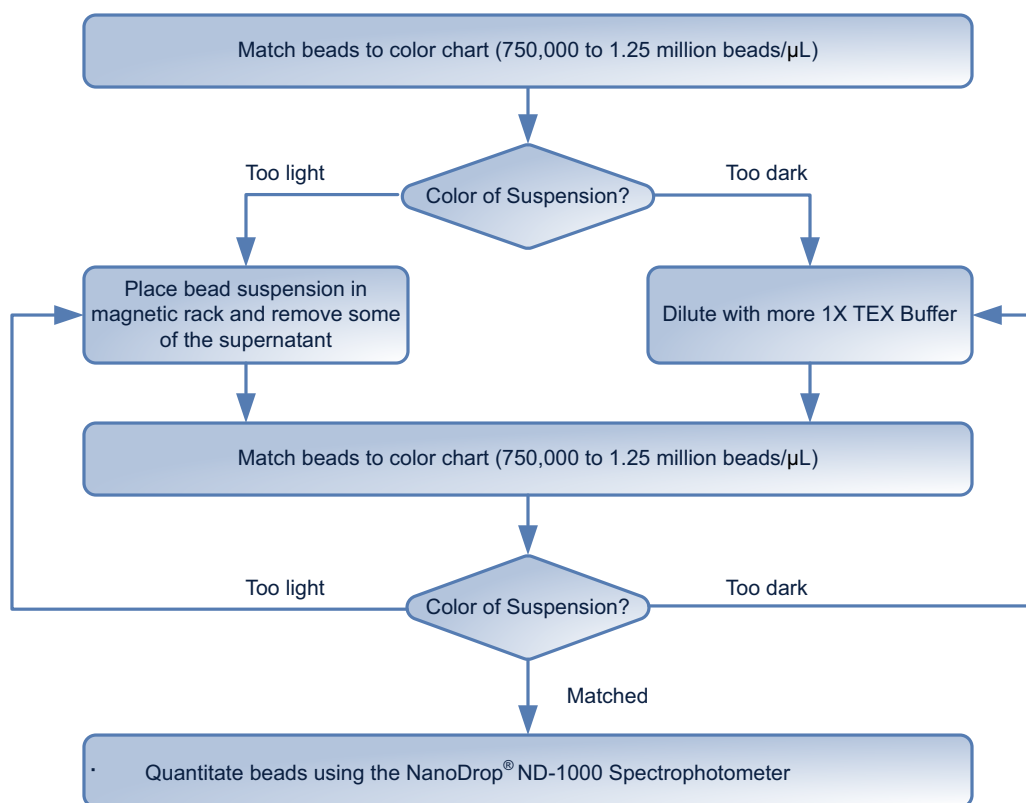


Figure 22 The SOLiD™ Bead Concentration Chart workflow.

5. When the bead concentration is within accurate range, quantitate the beads using the NanoDrop® ND-1000 Spectrophotometer. Take three readings, then average them. Calculate the bead concentration using the appropriate standard curve (for more details, see [“Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer”](#) on page 112).

STOPPING POINT. Store the templated beads at 4 °C in 1X TEX Buffer, or proceed to bead deposition and sequencing [refer to the *Applied Biosystems SOLiD™ 4 System Instrument Operation Guide* (4448379)].

4

Prepare Full-Scale Templated Beads

■ Prepare templated beads (full-scale)	46
■ Materials and equipment required (full-scale)	46
■ Workflow (full-scale)	46
■ Tips (full-scale)	47
■ Prepare the full-scale emulsion PCR (ePCR) reaction	49
■ Perform the emulsion break and bead wash (full-scale)	57
■ Enrich the full-scale templated beads	62
■ Modify the 3' ends (full-scale)	66

Prepare templated beads (full-scale)

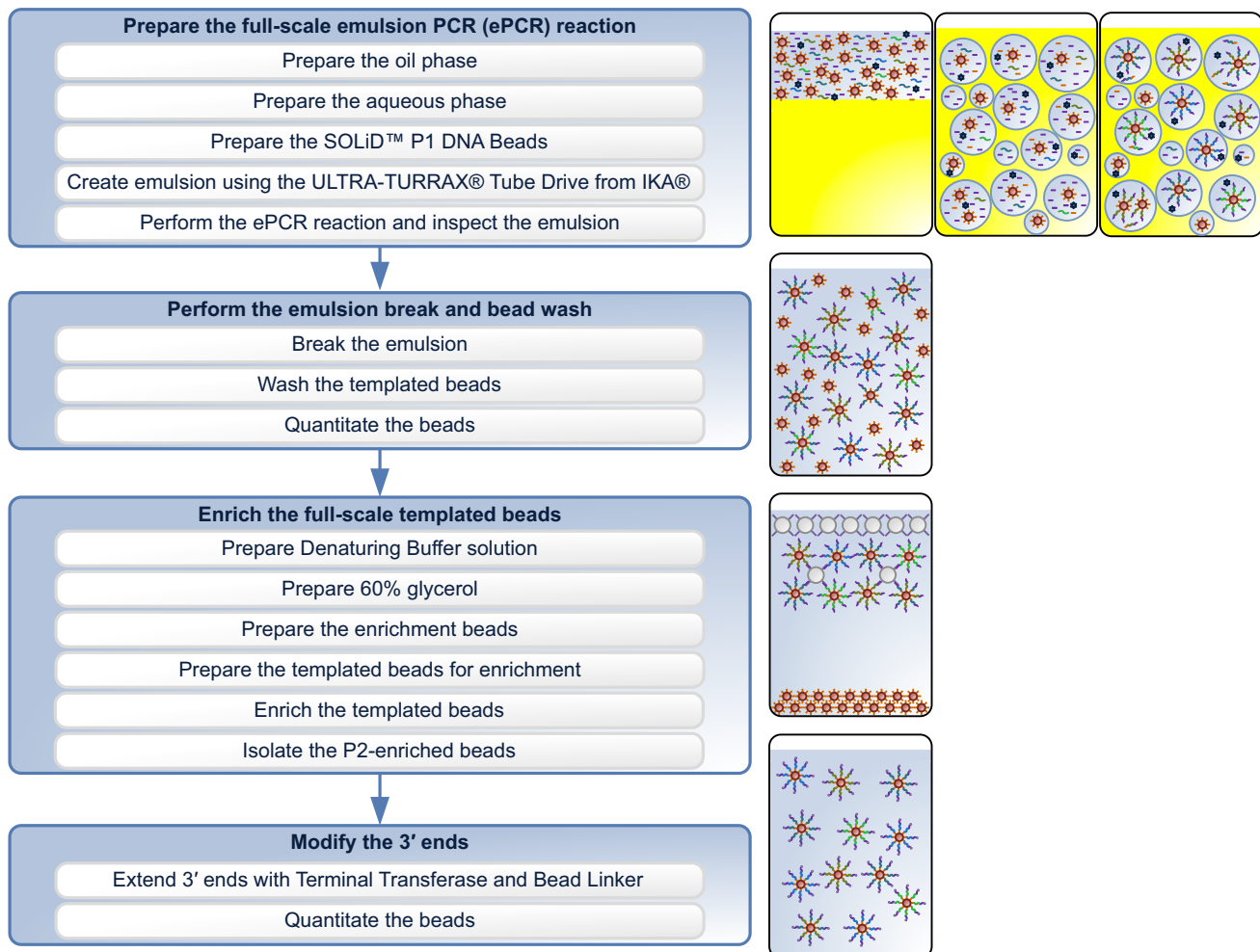
This chapter describes how to generate 150 to 300 million templated beads using the *full-scale* templated bead preparation method.

Materials and equipment required (full-scale)

See [Appendix A on page 95](#) for a list of equipment, kits, and consumables necessary for this procedure.

Workflow (full-scale)

See the descriptions of the steps below the workflow.



**Workflow overview
(full-scale)****Prepare the full-scale emulsion PCR (ePCR) reaction**

The oil phase and aqueous phase of the emulsion are prepared separately, then emulsified using the ULTRA-TURRAX® Tube Drive from IKA®. Each emulsion is seeded with 1.6 billion SOLiD™ P1 DNA Beads, then transferred into a single, 96-well plate for cycling. Different library template lengths require different numbers of cycles for thermal cycling.

Perform the emulsion break and bead wash (full-scale)

The emulsion break uses 2-butanol to purify emulsified templated beads from the oil phase following amplification. The beads are washed to remove residual 2-butanol, oil, and aqueous phase containing PCR reagents. The SOLiD™ Emulsion Collection Tray is placed over the 96-well plate, then the plate is centrifuged. Centrifuging the plate forces the emulsion from each well to a single reservoir. After centrifugation, 2-butanol is added to the reservoir. The broken emulsion is transferred to a 50-mL tube for further processing.

Enrich the full-scale templated beads

The templated bead enrichment procedure isolates beads with full-length extension products following ePCR. Beads with full-length extension products are isolated by oligo hybridization using the sequence of the P2 primer. Both monoclonal and polyclonal beads are enriched. The procedure is designed to enrich for templated beads derived from one full-scale ePCR reaction yielding 150 to 300 million templated beads.

Modify the 3' ends (full-scale)

The P2-enriched beads are extended with a Bead Linker by Terminal Transferase.

Tips (full-scale)

General

- Use syringes to accurately measure viscous reagents. Aspirate the volume very slowly from the reagent bottle to ensure that no air bubbles are trapped within the syringe. As is the best practice, draw some reagent into the syringe, dispense the entire reagent back to the reagent bottle, then draw the correct volume of reagent. Measure the volume to the point where the plunger contacts the side of the syringe.
- Perform all steps requiring 0.5-mL, 1.5-mL, and 2.0-mL tubes with Eppendorf LoBind Tubes. LoBind tubing from other vendors may have a chemical coating that has an adverse effect on bead deposition.
- Adjust microcentrifuge speeds and times according to the g-forces specified in the protocols. Applied Biosystems recommends the Eppendorf 5417R tabletop microcentrifuge.

**SOLiD™ P1 DNA
Beads**

- Do not freeze SOLiD™ P1 DNA Beads or templated beads. Store the SOLiD™ P1 DNA Beads at 4 °C in 1X TEX Buffer.

- If beads remain in the original tube after transfer, you can use a small additional volume of the appropriate buffer to recover the remaining beads. Do not exceed a total volume of 1.3 mL for a 1.5-mL LoBind Tube.

Covaris™ S2 System

- Applied Biosystems optimizes the procedures for the Covaris™ S2 System. The Covaris™ S2 System must be specially adapted to prepare beads for the Applied Biosystems SOLiD™ 4 System. Do not use the Covaris S1 sonicator or an unadapted Covaris™ S2 System for bead preparation. For more information, contact an Applied Biosystems SOLiD™ System applications specialist.
- To achieve optimal sonication by the Covaris™ S2 System, follow these guidelines:
 - Ensure that the Covaris™ S2 System is degassed.
 - Ensure that no bubbles are present in the Covaris™ S2 System.
 - Ensure that the instrument and tube are properly aligned for appropriate sonication of beads.
 - Use the appropriate adaptor with the Covaris™ S2 System, as shown in [Table 9](#). Place the tube collar at the indicator line of the adaptor.

Table 9 Tubes and adaptors for use with the Covaris™ S2 System

Sample volume	Size of LoBind Tube	Size of tube adaptor
≤ 200 µL	0.5-mL	0.65-mL
200 µL - 600 µL	1.5-mL	1.5-mL
600 µL - 1.2 mL	2.0-mL	1.5-mL

Prepare the full-scale emulsion PCR (ePCR) reaction

Prepare the oil phase (full-scale)

1. Use a *3-mL syringe* to dispense 1.8 mL of Emulsion Stabilizer 1 into the 50-mL conical tube.
2. Use a *1-mL syringe* to dispense 400 μL of Emulsion Stabilizer 2 very slowly into the 50-mL tube.
3. Pour the Emulsion Oil (approximately 37.8 mL) into the tube that has the Emulsion Stabilizer 1 and Emulsion Stabilizer 2 so that the final volume is 40 mL.
4. Cap the 50-mL tube, then vortex until all Emulsion Stabilizer 1 and Emulsion Stabilizer 2 are incorporated into the Emulsion oil.
5. Allow the mixture to degas for a minimum of 20 minutes while you prepare the aqueous phase (see [“Prepare the aqueous phase \(full-scale\)”](#)). To degas, place the mixture in a conical tube rack and slightly unscrew the conical tube cap.
6. Prime a 10-mL syringe by drawing in about 2 mL of oil phase then dispensing it back into the tube.
7. Take off the cap of a new SOLiD™ ePCR Tube. Use the primed 10-mL syringe to dispense 9 mL of oil phase into the SOLiD™ ePCR Tube, then cap the tube.

STOPPING POINT. The oil phase may be stored at 4 °C for up to 2 months. Before using the stored oil phase, thoroughly vortex and degas the solution for 20 minutes.

Prepare the aqueous phase (full-scale)

1. Dilute ePCR Primer 1 to prepare a 10- μM working stock solution. For each ePCR reaction, add 4 μL of ePCR Primer 1 to 36 μL of 1 \times Low TE buffer. Mix well.
2. Using only 1 \times Low TE Buffer and LoBind Tubes, prepare a dilution of the library template to a final concentration of 500 pM. Use [Table 10 on page 49](#) to convert the mass/volume concentration to molar concentration for each library (for calculation details, see [“Library Concentration Conversion” on page 117](#)). Dilute only enough template for the desired number of emulsions. If needed, perform a serial dilution of the library to accurately obtain the desired library concentration. For example, perform a 5 \times dilution from 50 nM to 10 nM, then perform a 20 \times dilution from 10 nM to 500 pM.

Table 10 Concentration conversions by library type

Library type	Average library length (bp)	Molar concentration (pM)	Mass/volume concentration
Fragment Library	~215	500	71 pg/ μL
2 x 50 bp Mate-Paired Library	~300	500	99 pg/ μL
2 x 25 bp Mate-Paired Library	~155	500	51 pg/ μL
Barcoded Fragment Library	~255	500	84 pg/ μL
Whole Transcriptome Library	~230	500	76 pg/ μL

Table 10 Concentration conversions by library type (*continued*)

Library type	Average library length (bp)	Molar concentration (pM)	Mass/volume concentration
Small RNA Library	~123	500	41 pg/μL
Barcoded SAGE Library	~130	500	43 pg/μL

⚠ **IMPORTANT!** Do not freeze-thaw dilutions of the library more than 3 to 4 times. Stock solutions and dilutions of libraries should be stored at – 20 °C at a concentration of 5 ng/μL or greater.

- Choose the appropriate library concentration, then prepare the aqueous phase by combining the following reagents in a Nalgene wide-mouth jar according to the table below (see [Table 11](#)). For workflow analysis, prepare the aqueous phase for library concentrations of 0.5 pM and 1.0 pM.

Table 11 Prepare the aqueous phase

		Library concentration		
Component	Final concentration [‡]	0.5 pM	1.0 pM	X pM
		Volume per reaction (μL) [§]		
10X PCR Buffer	1X	560	560	560
dNTP Mix (100 mM mix comprised of 25 mM each of dATP, dTTP, dCTP, dGTP)	14 mM (3.5 mM of each dNTP)	784	784	784
Magnesium Chloride (1 M)	25 mM	140	140	140
ePCR Primer 1 (10 μM working stock solution)	40 nM	22.4	22.4	22.4
ePCR Primer 2 (500 μM)	3 μM	33.6	33.6	33.6
Template (500 pM)	0.5 pM or 1.0 pM	5.6	11.2	X × 11.2
Nuclease-free Water	—	3294.4	3288.8	3300 – (X × 11.2)
AmpliTaq Gold® DNA Polymerase, UP (5 U/μL)	0.54 U/μL	600	600	600
Total	—	5440	5440	5440

[‡] The final concentration is based on a total volume of 5600 μL, which includes 5440 μL of liquid components and 160 μL of beads.

[§] Volumes below are for a single IKA®-based ePCR reaction to fill a 96-well plate.

- Keep the aqueous phase on ice until ready to use.

Prepare the SOLiD™ P1 DNA Beads (full-scale)

- Thoroughly vortex one tube of SOLiD™ P1 DNA Beads. Invert the tube at least once during vortexing to ensure that any beads stuck to the cap are washed down, then pulse-spin the tube.

2. Place the tube of beads in the magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
3. Resuspend the beads in 200 μ L of Bead Block Solution. Vortex the solution to ensure that all beads are suspended, then pulse-spin the tube.

⚠ **IMPORTANT!** Keep the Bead Block Solution at 4 °C until ready for use.

4. Sonicate the beads using the Bead Block Declump program on the Covaris™ S2 System (for program conditions, see [“Bead Block Declump” on page 127](#)), then pulse-spin the tube.
5. Place the tube in a magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
6. Resuspend the beads in 200 μ L of 1× TEX Buffer and vortex to ensure that all beads are suspended, then pulse-spin the beads.

Create the
emulsion with the
ULTRA-TURRAX®
Tube Drive from
IKA® (full-scale)

1. Place the SOLiD™ ePCR Tube containing 9 mL of oil phase onto the ULTRA-TURRAX® device, then twist the tube to lock it into position (see [Figure 23](#)).

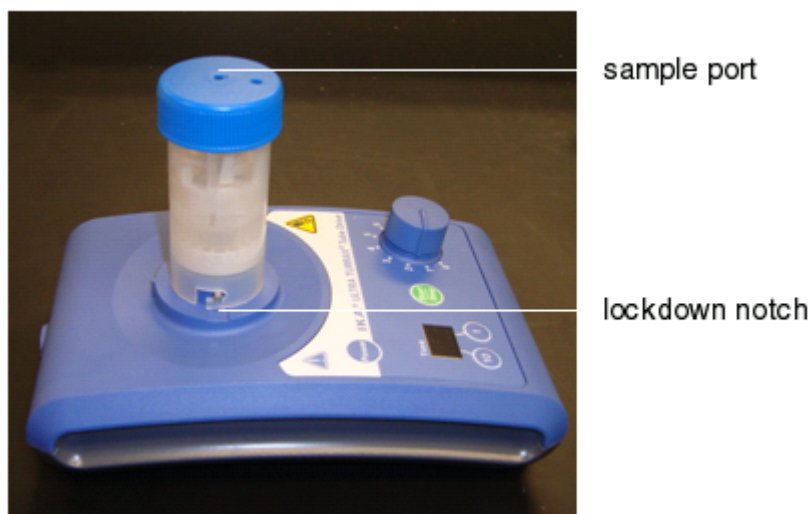


Figure 23 SOLiD™ ePCR Tube on the ULTRA-TURRAX® Tube Drive from IKA®.

2. Sonicate the SOLiD™ P1 DNA beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)), then pulse-spin the beads.
3. Immediately add 160 μ L of the SOLiD™ P1 DNA Beads to the aqueous phase, then mix by gently swirling the bottle to ensure that the beads are uniformly dispersed (see [Figure 24](#)).



Figure 24 SOLiD™ P1 DNA Beads mixed in aqueous phase.

4. Verify that the Xstream pipettor is set up for full-scale emulsions (see [Figure 25](#)):
 - Dial Setting: **Pip**
 - Speed (aspirate UP): **scale 5** (mid-range)
 - Speed (dispense DOWN): **scale 1** (slowest)
 - Total volume: **5.60 mL**

If necessary, reprogram the Xstream pipettor (see [“Program the Eppendorf Repeater® Xstream Pipettor” on page 106](#)).



Figure 25 Xstream pipettor settings, with the dial setting set to Pip.

5. Attach a 10-mL Combitip Plus tip onto the Xstream pipettor.
6. Fill the 10-mL Combitip Plus tip with the entire 5.60 mL of aqueous phase and bead mixture with the Xstream pipettor (see [Figure 26 on page 53](#)).



Figure 26 Filling the 10-mL Combipip Plus tip with the aqueous phase and bead mixture using the Xstream pipettor.

7. Verify that the time on the ULTRA-TURRAX[®] Tube Drive from IKA[®] is set to **5** minutes, then press the **Start** button.
8. Wait for the instrument's fly wheel to engage and to reach proper speed, then gently place the Combipip Plus tip into the center sample loading hole in the ULTRA-TURRAX[®] cap (see [Figure 27 on page 54](#)).



Figure 27 Correct placement of Combitip Plus into sample port in SOLiD™ ePCR Tube cap.

9. Dispense the aqueous phase and bead mixture into the spinning oil phase. When the entire volume is dispensed, press the center blue button *twice* on the pipettor to empty all contents from the Combitip Plus tip.
10. Remove a 5-mL Combitip Plus tip from its packaging, then cut off its end at the bevel with a razor blade (see [Figure 28 on page 55](#)).

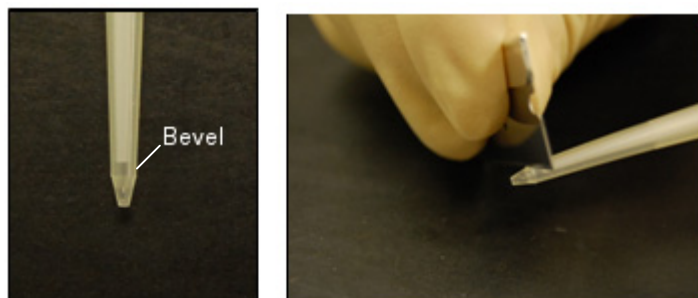


Figure 28 Cutting the Combipip Plus tip for emulsion dispersion.

11. Attach the cut Combipip Plus tip onto an Eppendorf Repeater® Plus Pipette.
12. Gently dispense 150 μ L of emulsion into each well of a 96-well PCR plate, then seal the plate with clear adhesive film (see Figure 29).



Figure 29 Emulsion transferred to a 96-well plate.

Perform the ePCR reaction and inspect the emulsion (full-scale)

1. Set up the ePCR conditions on the GeneAmp® PCR System 9700:
 - ePCR thermal cycling program:

Stage	Step	Temp (°C)	Time
Holding	Denature	95	5 min
40 cycles [‡] or 60 cycles [§]	Denature	93	15 sec
	Anneal	62	30 sec
	Extend	72	75 sec
Holding	Final extension	72	7 min
Holding	—	4	∞

[‡] Set 40 cycles: Fragment library or 2 \times 25 bp mate-paired library.

[§] Set 60 cycles: 2 \times 50 bp mate-paired library.

- Ramp speed: **9600**
- Reaction volume: **50 μ L**

2. Place the 96-well plate in a GeneAmp® PCR System 9700, then start the run.
3. After the ePCR Program finishes, inspect the bottom of the reaction plate for beads that have fallen out of the emulsion. Beads appear as amber-colored specks at the bottom of a well. A small number of beads may fall out of emulsion and appear as small brown flecks at the bottom of a well. Applied Biosystems does not recommend any further processing of emulsions that have more than 3 wells of broken emulsion, where aqueous phase appears at the bottom of a well (see [Figure 30](#)).

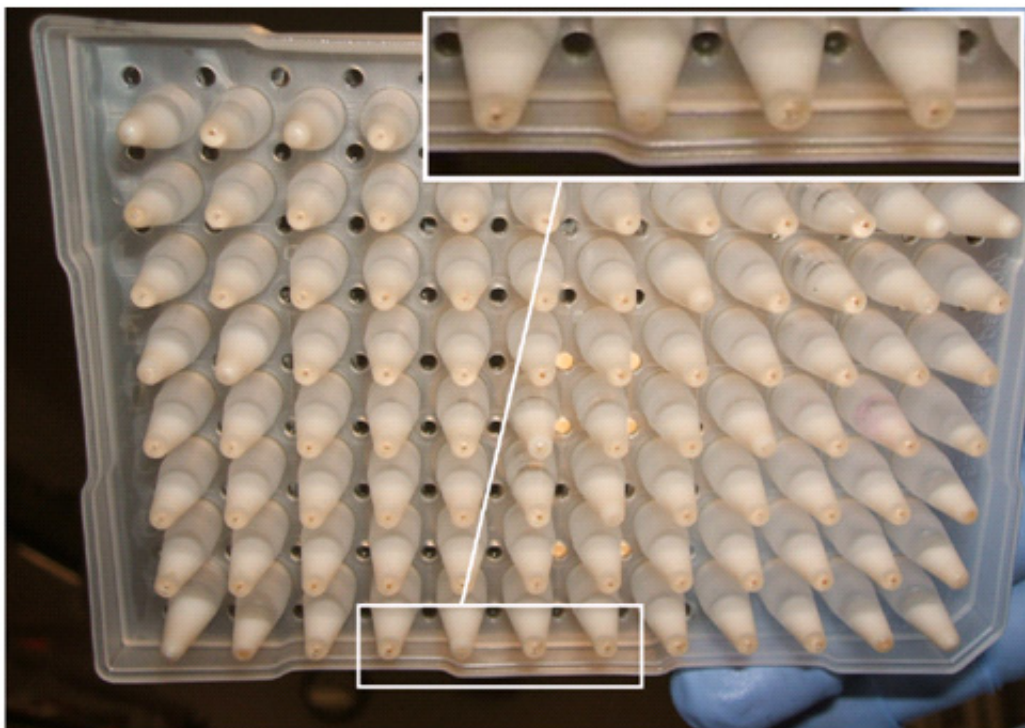


Figure 30 Broken emulsions.

STOPPING POINT. Store the 96-well plate at 4 °C, or proceed to [“Perform the emulsion break and bead wash \(full-scale\)” on page 57](#).

Perform the emulsion break and bead wash (full-scale)

Break the emulsion (full-scale)

1. Place the SOLiD™ Emulsion Collection Tray on top of the ePCR 96-well plate (see [Figure 31](#)).

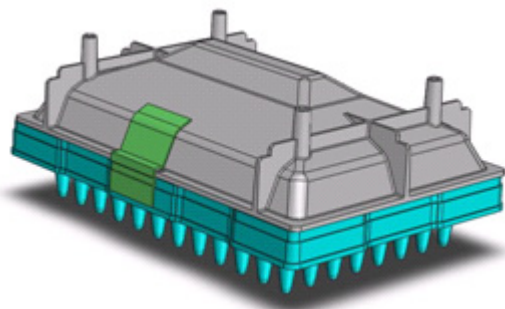


Figure 31 The SOLiD™ Emulsion Collection Tray taped to a 96-well reaction plate.

2. Seal the pieces together with tape on all four sides. Immediately prior to centrifugation, flip the entire apparatus so that the 96-well plate is upside-down over the collection tray (see [Figure 32](#)).

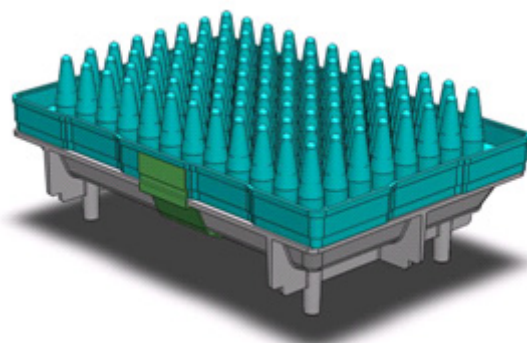


Figure 32 The inverted plate assembly.

3. Centrifuge the inverted plate and reservoir for 2 minutes at $550 \times g$ according to the recommended centrifuge settings (see [Table 12](#)).

Table 12 Recommended centrifuge settings

Adjustable parameter	Recommend setting
Acceleration (independent)	High
Deceleration (independent)	Low
Acceleration/Deceleration (single setting)	Mid

4. After centrifugation is complete, remove the plate assembly from the centrifuge and place on the lab bench. Hold the assembly steady, then gently remove the tape and 96-well plate from the collection tray.



Note: Ensure that the centrifuge is working properly and maintained regularly. Use anti-slip pads in the centrifuge carriers whenever possible.

5. In a fume hood, add 10 mL of 2-butanol to the collection tray using a serological pipette.
6. Pipet the emulsion up and down until the solution is homogeneous.
7. Transfer all the emulsion and 2-butanol to a 50-mL conical tube.
8. Rinse the reservoir with an additional 6 mL of 2-butanol to ensure that all residual beads are recovered.
9. Cap the tube, then vortex to mix the solution.
10. Centrifuge the tube at $2000 \times g$ for 5 minutes.
11. Gently decant the 2-butanol-oil phase into a waste bottle. With the tube inverted, place the tube onto paper towels to drain residual 2-butanol-oil.
12. Wait 5 minutes to ensure that all the oil is removed.



IMPORTANT! If the pellet begins to slide out, stop decanting, then remove the 2-butanol using a pipette.

Wash the templated beads (full-scale)

1. Place the 50-mL tube upright in a rack, then add 600 μ L of 1 \times Bead Wash Buffer. Let the pellet soak in 1 \times Bead Wash Buffer for 2 minutes.
2. Resuspend the pellet by gently pipetting up and down, then transfer the beads from the 50-mL tube to a 1.5-mL LoBind Tube.
3. Rinse the *bottom* of the 50-mL tube with an additional 600 μ L of 1 \times Bead Wash Buffer, then transfer the wash to the 1.5-mL LoBind Tube.
4. Vortex the 1.5-mL LoBind Tube, then centrifuge the tube at $21,000 \times g$ (minimum $14,000 \times g$) for 1 minute.
5. Remove the top oil phase with a pipette. Remove as much of the oil at the meniscus as possible.
6. With a new pipette tip, carefully remove and discard the supernatant.
7. Resuspend the pellet by adding 150 μ L of 1 \times Bead Wash Buffer to the tube, then vortex the tube. Pulse-spin the tube, then transfer the mixture to a new 1.5-mL LoBind Tube.
8. Rinse the *bottom* of the original tube with an additional 150 μ L of 1 \times Bead Wash Buffer, then transfer the wash to the new tube.
9. Add 1 mL of 1 \times Bead Wash Buffer to the new tube, then vortex the tube.
10. Centrifuge the tube at $21,000 \times g$ (minimum $14,000 \times g$) for 1 minute, then remove and discard the supernatant.
11. Resuspend the beads in 200 μ L of 1 \times TEX Buffer.
12. Place the tube in a magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
13. Resuspend the beads in 200 μ L of 1 \times TEX Buffer.

STOPPING POINT. Store the templated beads at 4 °C in 1 \times TEX Buffer, or proceed to [“Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer \(full-scale\)”](#) or to [“Quantitate the beads using a hemocytometer”](#) on page 115.

Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer (full-scale)

1. If necessary, generate a standard curve (see [“Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer”](#) on page 112).
2. Sonicate the beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1”](#) on page 127), then pulse-spin the beads.

3. Make a 1-mL dilution of beads in 1× TEX Buffer (1:10 dilution recommended) in a 1.5-mL LoBind Tube.
4. Use the SOLiD™ Bead Concentration Chart (Applied Biosystems PN 4415131) to estimate the bead concentration of the beads (see Figure 33 for a picture of the chart; for best results use the official chart).

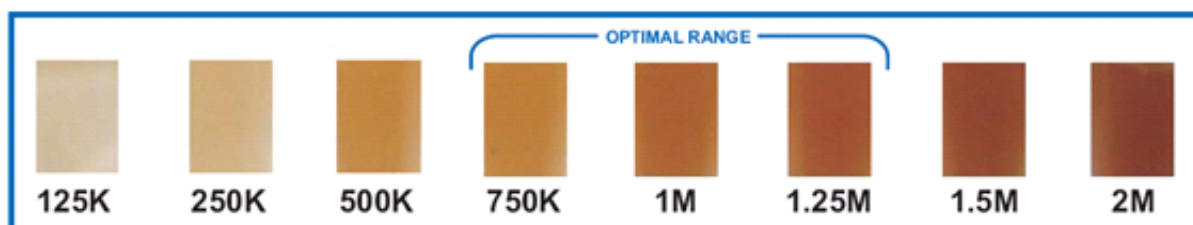


Figure 33 The SOLiD™ Bead Concentration Chart (facsimile). For best results, use the SOLiD™ Bead Concentration Chart (PN 4415131), supplied separately.

5. Adjust the volume of beads so that the color of the bead solution matches a color in the optimal range (750,000 beads/μL to 1.25 million beads/μL; see Figure 34 for the workflow).

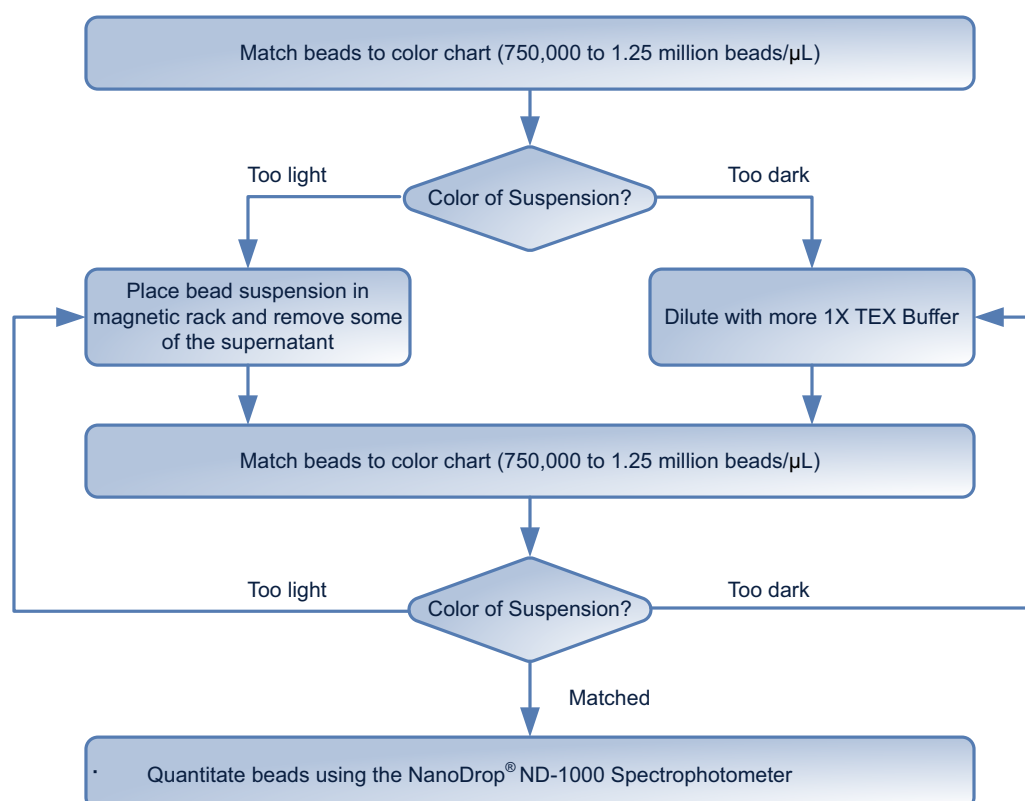


Figure 34 The SOLiD™ Bead Concentration Chart workflow.

6. When the bead concentration is within accurate range, quantitate the beads using the NanoDrop® ND-1000 Spectrophotometer. Take three readings, then average them. Calculate the bead concentration using the appropriate standard curve (for more details, see [“Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer”](#) on page 112).
7. Combine all diluted and undiluted beads.

STOPPING POINT. Store the beads at 4 °C in 1X TEX Buffer, or proceed to [“Enrich the full-scale templated beads”](#) on page 62.

Enrich the full-scale templated beads

Prepare the Denaturing Buffer solution (full-scale)

1. For each ePCR reaction, transfer 1.8 mL of Denaturing Buffer to a 15-mL conical tube.
 2. Add 200 μ L of Denaturant to the 1.8 mL of Denaturing Buffer, then cap the tube and vortex.
- !** **IMPORTANT!** Create a new batch of the prepared Denaturing Buffer solution each week.

Prepare 60% glycerol (full-scale)

1. With a 10-mL syringe, add 4 mL of Nuclease-free Water to a 15-mL conical tube.
 2. With a 3-mL syringe, add 6 mL glycerol to the Nuclease-free Water by dispensing 3 mL glycerol twice with the syringe. Fill and dispense the glycerol slowly to ensure that the total volume of glycerol is dispensed.
 3. Cap the tube, then vortex to mix the solution well.
- !** **IMPORTANT!** Prepare the 60% glycerol fresh weekly.

Prepare the enrichment beads (full-scale)

1. Vortex the enrichment beads and immediately transfer 650 μ L of the enrichment beads to a 1.5-mL LoBind Tube.
2. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
3. Resuspend the enrichment beads in 900 μ L of 1 \times Bind & Wash Buffer.
4. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
5. Repeat steps 3 and 4.
6. Resuspend the enrichment beads in 350 μ L of 1 \times Bind & Wash Buffer.
7. Add 3.5 μ L of 1 mM Enrichment Oligo to the tube of enrichment beads, then vortex and pulse-spin the tube.
8. Rotate the tube at room temperature for 30 minutes.
9. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove the supernatant.
10. Resuspend the enrichment beads in 900 μ L of 1 \times TEX Buffer.

11. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
12. Repeat steps 10 and 11.
13. Resuspend the enrichment beads in 150 μL of 1 \times Low Salt Binding Buffer.

STOPPING POINT. Store the prepared enrichment beads at 4 °C in 1 \times Low Salt Binding Buffer, or proceed to [“Prepare the templated beads for enrichment \(full-scale\)”](#). Prepared enrichment beads should be used within one week of preparation.

**Prepare the
templated beads
for enrichment (full-
scale)**

1. Place the tube in the magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
2. Resuspend the templated beads in 300 μL of prepared Denaturing Buffer solution, then let the suspension stand for 1 minute.
3. Place the tube in the magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
4. Repeat steps 2 and 3 twice.
5. Resuspend the templated beads in 300 μL of 1 \times TEX Buffer.
6. Place the tube in the magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
7. Repeat steps 5 and 6 twice.
8. Resuspend the templated beads in 150 μL of 1 \times TEX Buffer, then transfer the templated bead suspension to a new 0.5-mL LoBind Tube.
9. Sonicate the beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)).

Enrich the templated beads (full-scale)

1. Transfer the prepared enrichment beads to the tube of templated beads, then vortex and pulse-spin the bead mixture.
2. Sonicate the enrichment-templated bead mixture using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 3” on page 127](#)), then pulse-spin the beads.
3. Incubate the bead mixture at 61 °C for 15 minutes. During the incubation, vortex and pulse-spin the bead mixture every 5 minutes including at the end of the incubation.
4. Immediately cool the beads on ice for 2 minutes.
5. Add 600 µL of *freshly prepared* 60% glycerol to a new 1.5-mL LoBind Tube.
6. Gently pipet the bead mixture up and down the beads to mix, then load the entire volume *carefully* on top of the 60% glycerol solution. Do *not* vortex the tube.
7. Centrifuge the tube for 3 minutes at $21,000 \times g$ (minimum $14,000 \times g$).
8. Add 1 mL of 1× TEX Buffer to a new 2.0-mL LoBind Tube.
9. Transfer the top layer of beads into the tube with 1× TEX Buffer. Aspirate as little glycerol as possible to collect all of the beads at the top layer without touching the un-templated beads at the bottom of the tube. When you dispense the top layer of beads into the 1× TEX Buffer, dispense the beads into the bottom of the tube. Aspirate a small amount of 1× TEX buffer to clean the pipette tip.
10. Top off the tube with additional 1× TEX Buffer to the 2.0-mL mark, then vortex.
11. Centrifuge the tube for 1 minute at $21,000 \times g$ (minimum $14,000 \times g$).



Note: Verify the beads are pelleted in case excess glycerol carried over to the 1× TEX Buffer creates a matrix that impedes pelleting of beads.

Proceed according to the table below ([Table 13](#)).

Table 13 Steps for pelleted or unpelleted beads


If the beads are...	Then perform steps...
Pelleted	12 and 13
Not pelleted	14 to 16

12. Remove the supernatant. Add 400 µL of 1× TEX Buffer to the tube of beads and vortex.
13. Proceed to [“Isolate the P2-enriched beads \(full-scale\)” on page 65](#).
14. Transfer half of the tube volume to a new 2.0-mL LoBind Tube, then add an additional 500 µL of 1× TEX Buffer to each tube. Vortex each tube.

15. Centrifuge the tubes for 1 minute at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
16. Add 200 μL of 1 \times TEX Buffer to each tube, resuspend the beads, then pool the beads into one tube.

**Isolate the
P2-enriched beads
(full-scale)**

1. Centrifuge the tube for 1 minute at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.

 **IMPORTANT!** *Never* magnet the P2-enriched beads before adding prepared Denaturing Buffer solution to the beads. If you do, the templated beads linked to the enrichment beads are lost when the supernatant is removed.
2. Resuspend the pellet with 400 μL of prepared Denaturing Buffer solution, then let stand for 1 minute.
3. Place the tube in a magnetic rack for at least 1 minute until the supernatant is pure white or clear, then remove and discard the supernatant.
4. Repeat steps 2 and 3 until the supernatant is clear (all white enrichment beads have been removed).
5. Resuspend the beads in 400 μL of 1 \times TEX Buffer.
6. Place the tube in a magnetic rack for at least 1 minute until the supernatant is clear. Remove and discard the supernatant.
7. Repeat steps 5 and 6.
8. Resuspend the beads in 200 μL 1 \times TEX Buffer. Vortex, pulse-spin, then transfer the bead solution to a 1.5-mL LoBind Tube.
9. Rinse the 2.0-mL tube with 200 μL 1 \times TEX Buffer and transfer the rinse to the 1.5-mL LoBind Tube.
10. Sonicate the beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)). Pulse-spin the beads.
11. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
12. Resuspend the beads in 400 μL of 1 \times TEX Buffer.
13. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
14. If the supernatant appears cloudy due to residual enrichment beads, repeat steps 12 and 13 until the supernatant is clear.

15. Resuspend the beads in 400 μL of 1 \times TEX Buffer.

STOPPING POINT. Store the prepared enrichment beads at 4 $^{\circ}\text{C}$ in 1 \times TEX Buffer, or proceed to “[Modify the 3' ends \(full-scale\)](#)”.

Modify the 3' ends (full-scale)

Extend the 3' ends with Terminal Transferase and Bead Linker (full- scale)

1. If the P2-enriched beads have been stored overnight or longer, sonicate the beads using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see “[Covalent Declump 3](#)” on page 127). Pulse-spin the beads.
2. For each ePCR reaction, prepare 500 μL of 1 \times Terminal Transferase Reaction Buffer according to [Table 14](#):

Table 14 Prepare 1 \times Terminal Transferase Reaction Buffer

Component	Volume per reaction (μL)
10 \times Terminal Transferase Buffer	55
10 \times Cobalt Chloride	55
Nuclease-free Water	390
Total	500



Note: The 1 \times Terminal Transferase Reaction Buffer should be clear. If the solution becomes colored, discard then prepare fresh buffer using a new lot of material.

3. Add 1 μL of 50 mM Bead Linker to 49 μL of 1 \times Low TE Buffer to prepare a 1 mM Bead Linker solution.
4. Place the tube of P2-enriched beads in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
5. Resuspend the beads in 100 μL of 1 \times Terminal Transferase Reaction Buffer, then transfer the beads to a new 1.5-mL LoBind Tube.
6. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
7. Resuspend the beads in 100 μL of 1 \times Terminal Transferase Reaction Buffer.
8. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
9. Resuspend the beads in 178 μL of 1 \times Terminal Transferase Reaction Buffer.
10. Add 20 μL of 1 mM Bead Linker solution.

11. Sonicate the beads using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 3” on page 127](#)). Pulse-spin the beads.
12. Add 2 μL of Terminal Transferase (20 U/ μL) and vortex. Pulse-spin the beads.
13. Place the tube on a rotator and rotate for 2 hours at 37 °C.
14. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
15. Resuspend the beads in 400 μL of 1× TEX Buffer.
16. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
17. Resuspend the beads in 400 μL of 1× TEX Buffer.

STOPPING POINT. Store the templated beads at 4 °C in 1× TEX Buffer, or proceed to [“Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer \(full-scale\)”](#) or [“Quantitate the beads using a hemocytometer” on page 115](#).

Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer (full-scale)

1. If necessary, generate a standard curve (see [“Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer” on page 112](#)).
2. Sonicate the beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)), then pulse-spin the beads.
3. Use the SOLiD™ Bead Concentration Chart (Applied Biosystems PN 4415131) to estimate the bead concentration of the beads (see [Figure 35](#) for a picture of the chart; for best results use the official chart).

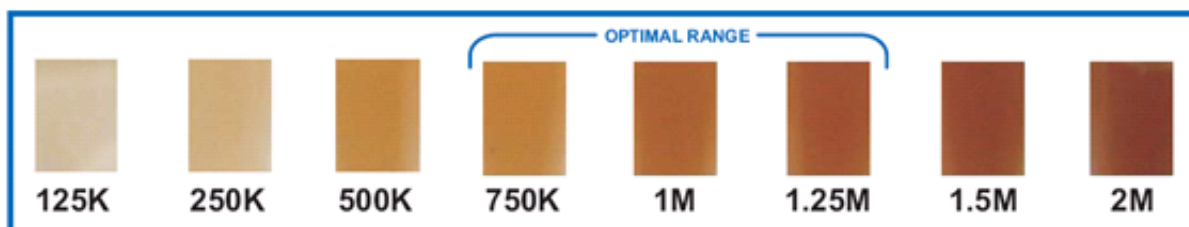


Figure 35 The SOLiD™ Bead Concentration Chart (facsimile). For best results, use the SOLiD™ Bead Concentration Chart (PN 4415131), supplied separately.

4. Adjust the volume of beads so that the color of the bead solution matches a color in the optimal range (750,000 beads/ μL to 1.25 million beads/ μL ; see [Figure 36 on page 68](#) for the workflow).

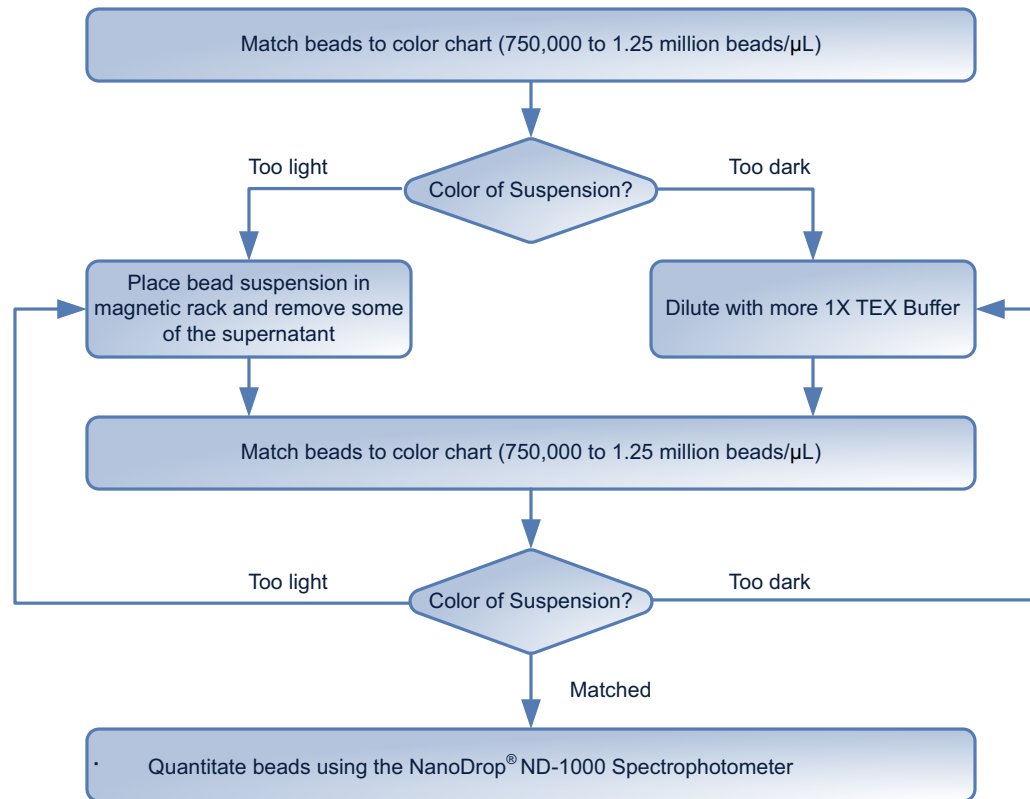


Figure 36 The SOLiD™ Bead Concentration Chart workflow.

5. When the bead concentration is within accurate range, quantitate the beads using the NanoDrop® ND-1000 Spectrophotometer. Take 3 readings, then average them. Calculate the bead concentration using the appropriate standard curve (for more details, see [“Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer”](#) on page 112).

STOPPING POINT. Store the templated beads at 4 °C in 1X TEX Buffer, or proceed to bead deposition and sequencing [refer to the *Applied Biosystems SOLiD™ 4 System Instrument Operation Guide* (PN 4448379)].

5

Macro-Scale Templated Beads (4 and 8 ePCR Reactions)

- Prepare templated beads (macro-scale) 70
- Materials and equipment required (macro-scale) 70
- Workflow (macro-scale) 71
- Tips (macro-scale) 72
- Macro-scale emulsion instructions 74

Prepare templated beads (macro-scale)

This chapter describes preparation steps and the workflow overview for generating 600 million to 2.4 billion templated beads using the *macro*-scale templated bead preparation method.

After completing the instructions in this chapter, you must complete the instructions in either [Chapter 6](#), for 4 ePCR reactions, or [Chapter 7](#), for 8 ePCR reactions:

Macro-Scale (4 ePCR reactions)

The emulsion PCR protocol to prepare macro-scale (4 ePCR reactions) templated beads is described in [Chapter 6](#), “[Prepare Macro-Scale Templated Beads \(4 ePCR Reactions\)](#)” on page 75.

Macro-Scale (8 ePCR reactions)

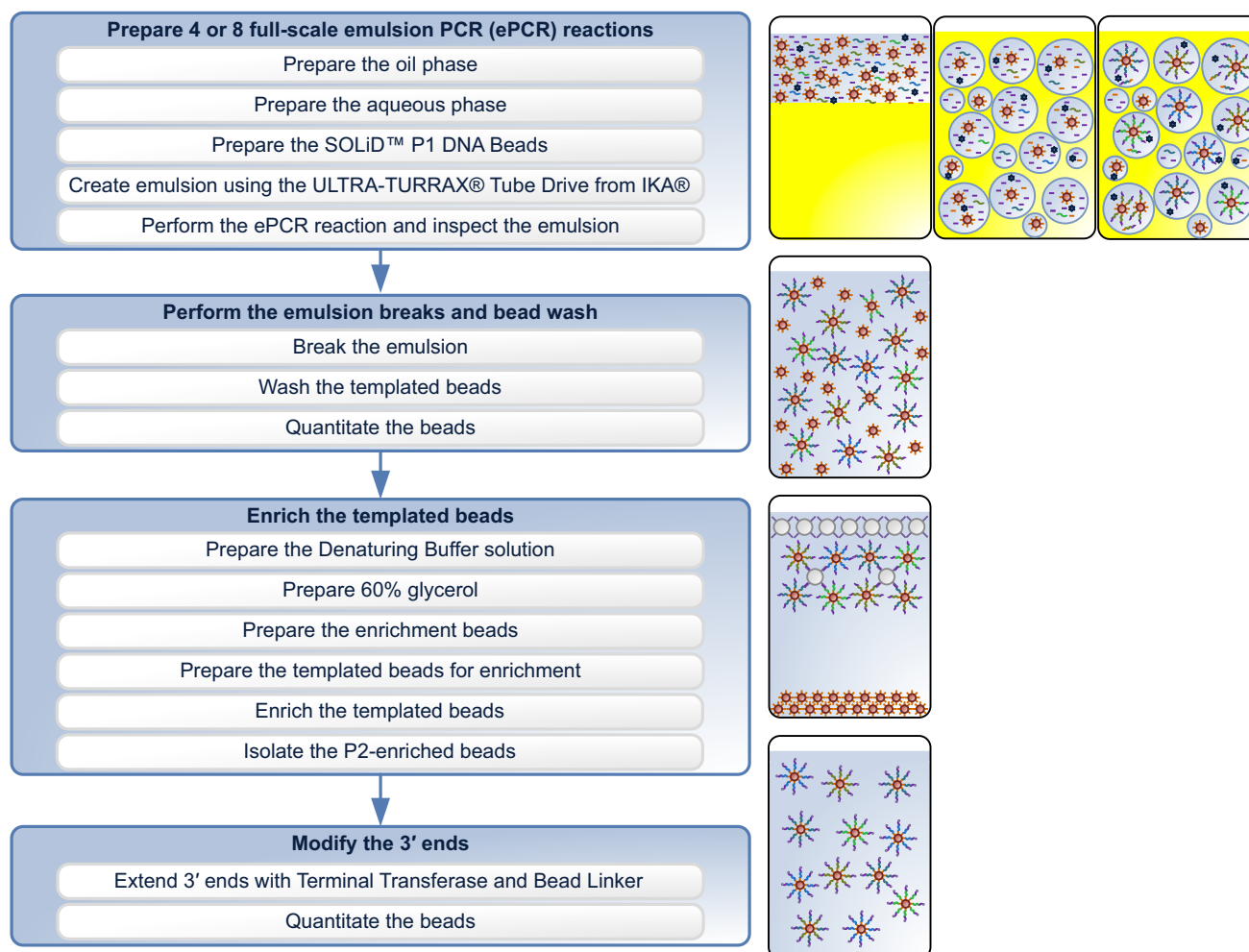
The emulsion PCR protocol to prepare macro-scale (8 ePCR reactions) templated beads is described in [Chapter 7](#), “[Prepare Macro-Scale Templated Beads \(8 ePCR Reactions\)](#)” on page 85.

Materials and equipment required (macro-scale)

See [Appendix A on page 95](#) for a list of equipment, kits, and consumables necessary for this procedure.

Workflow (macro-scale)

See the overview descriptions of the steps below the workflow.



Workflow overview (macro-scale)

Prepare 4 or 8 full-scale emulsion PCR (ePCR) reactions

The oil phase and aqueous phase of the emulsion are prepared separately, then emulsified using the ULTRA-TURRAX® Tube Drive from IKA®. Each emulsion is seeded with 1.6 billion SOLiD™ P1 DNA Beads, then transferred into a single, 96-well plate for cycling. Depending on the output required for your experiment, perform 4 or 8 ePCR reactions. Different library template lengths require different numbers of cycles for thermal cycling.

Perform the emulsion break and bead wash (macro-scale)

The emulsion break uses 2-butanol to purify emulsified templated beads from the oil phase following amplification. The beads are washed to remove any residual 2-butanol, oil, and aqueous phase containing PCR reagents. The SOLiD™ Emulsion Collection Tray is placed over the 96-well plate, then the plate is centrifuged. Centrifuging the plate forces the emulsion from each well to a single reservoir. After centrifugation, 2-butanol is added to the reservoir. The broken emulsion is transferred to a 50-mL tube for further processing.

Enrich the templated beads (macro-scale)

The templated bead enrichment procedure isolates beads with full-length extension products following ePCR. Beads with full-length extension products are isolated by oligo hybridization using the sequence of the P2 primer. Both monoclonal and polyclonal beads are enriched. The procedure is designed to enrich the templated beads derived from four or eight ePCR reactions containing 1.6 billion SOLiD™ P1 DNA Beads each (6.4 billion SOLiD™ P1 DNA Beads for four ePCR reactions or 12.8 billion SOLiD™ P1 DNA Beads for eight ePCR reactions).

Modify the 3' ends (macro-scale)

The P2-enriched beads are extended with a Bead Linker by Terminal Transferase.

Tips (macro-scale)**General**

- Use syringes to accurately measure viscous reagents. Aspirate the volume very slowly from the reagent bottle to ensure that no air bubbles are trapped within the syringe. As is the best practice, draw some reagent into the syringe, dispense the entire reagent back to the reagent bottle, then draw the correct volume of reagent. Measure the volume to the point where the plunger contacts the side of the syringe.
- Perform all steps requiring 0.5-mL, 1.5-mL, and 2.0-mL tubes with Eppendorf LoBind Tubes. LoBind tubing from other vendors may have a chemical coating that has an adverse effect on bead deposition.
- Adjust microcentrifuge speeds and times according to the g-forces specified in the protocols. Applied Biosystems recommends the Eppendorf 5417R tabletop microcentrifuge.

SOLiD™ P1 DNA Beads

- Do not freeze SOLiD™ P1 DNA Beads or templated beads. Store the SOLiD™ P1 DNA Beads at 4 °C in 1× TEX Buffer.
- If beads remain in the original tube after transfer, you can use a small additional volume of the appropriate buffer to recover the remaining beads. Do not exceed a total volume of 1.3 mL for a 1.5-mL LoBind Tube.

Covaris™ S2 System

- Applied Biosystems optimizes the procedures for the Covaris™ S2 System. The Covaris™ S2 System must be specially adapted to prepare beads for the Applied Biosystems SOLiD™ 4 System. Do not use the Covaris S1 sonicator or an unadapted Covaris™ S2 System for bead preparation. For more information, contact an Applied Biosystems SOLiD™ System applications specialist.
- To achieve optimal sonication by the Covaris™ S2 System, follow these guidelines:
 - Ensure that the Covaris™ S2 System is degassed.
 - Ensure that no bubbles are present in the Covaris™ S2 System.
 - Ensure that the instrument and tube are properly aligned for appropriate sonication of beads.
 - Use the appropriate adaptor with the Covaris™ S2 System, as shown in [Table 15](#). Place the tube collar at the indicator line of the adaptor.

Table 15 Tubes and adaptors for use with the Covaris™ S2 System™

Sample volume	Size of LoBind Tube	Size of tube adaptor
≤ 200 µL	0.5-mL	0.65-mL
200 µL - 600 µL	1.5-mL	1.5-mL
600 µL - 1.2 mL	2.0-mL	1.5-mL

Macro-scale emulsion instructions

After completing the instructions in this chapter, you must complete the instructions in either [Chapter 6, Prepare Macro-Scale Templated Beads \(4 ePCR Reactions\)](#) or [Chapter 7, Prepare Macro-Scale Templated Beads \(8 ePCR Reactions\)](#).

Macro-Scale (4 ePCR reactions)

The emulsion PCR protocol to prepare macro-scale (4 ePCR reactions) templated beads is described in [Chapter 6, “Prepare Macro-Scale Templated Beads \(4 ePCR Reactions\)”](#) on page 75.

Macro-Scale (8 ePCR reactions)

The emulsion PCR protocol to prepare macro-scale (8 ePCR reactions) templated beads is described in [Chapter 7, “Prepare Macro-Scale Templated Beads \(8 ePCR Reactions\)”](#) on page 85.

6

Prepare Macro-Scale Templated Beads (4 ePCR Reactions)

- Prepare the templated beads (macro-scale: 4 ePCR reactions). 76
- Prepare 4 full-scale emulsion PCR (ePCR) reactions. 76
- Perform the emulsion breaks and bead wash (macro-scale: 4 ePCR reactions). . 76
- Enrich the templated beads (macro-scale: 4 ePCR reactions) 76
- Modify the 3' ends (macro-scale: 4 ePCR reactions) 82

Prepare the templated beads (macro-scale: 4 ePCR reactions)

These instructions describe how to generate 600 million to 1.2 billion templated beads using the *macro-scale* (4 ePCR reactions) templated bead preparation method.

Note: Do not begin these instructions unless you have completed the instructions in [Chapter 5, Macro-Scale Templated Beads \(4 and 8 ePCR Reactions\)](#). [Chapter 5](#) also describes the workflow overview.

To instead prepare macro-scale (8 ePCR reactions), use [Chapter 7, “Prepare Macro-Scale Templated Beads \(8 ePCR Reactions\)”](#) on page 85.

Prepare 4 full-scale emulsion PCR (ePCR) reactions

Prepare 4 emulsion PCR reactions as described in “[Prepare the full-scale emulsion PCR \(ePCR\) reaction](#)” in [Chapter 4, Prepare Full-Scale Templated Beads](#). Four ePCR reactions provide an adequate bead yield for two wells of a 4-well slide for sequencing. Store each 96-well plate at 4 °C or proceed to “[Perform the emulsion breaks and bead wash \(macro-scale: 4 ePCR reactions\)](#)”.


Perform the emulsion breaks and bead wash (macro-scale: 4 ePCR reactions)

Perform the emulsion break and bead wash procedure on each of the 4 emulsion PCR reactions, as described in “[Perform the emulsion break and bead wash \(full-scale\)](#)” in [Chapter 4, Prepare Full-Scale Templated Beads](#). Store each tube of beads at 4 °C or proceed to “[Enrich the templated beads \(macro-scale: 4 ePCR reactions\)](#)”.

Enrich the templated beads (macro-scale: 4 ePCR reactions)

Prepare the Denaturing Buffer solution


1. For each enrichment (4 plates to be combined), transfer 5.4 mL of Denaturing Buffer to a 15-mL conical tube.
2. Add 600 µL of Denaturant to the 5.4 mL of Denaturing Buffer, then cap the tube and vortex.

 **IMPORTANT!** Create a new batch of the prepared Denaturing Buffer solution each week.

Prepare 60% glycerol

1. With a 10-mL syringe, add 4 mL of Nuclease-free Water to a 15-mL conical tube.

2. With a 3-mL syringe, add 6 mL of glycerol to the Nuclease-free Water by dispensing 3 mL of glycerol twice with the syringe. Fill and dispense the glycerol slowly to ensure that the total volume of glycerol is dispensed.
3. Cap the tube, then vortex to mix the solution well.

 **IMPORTANT!** Prepare the 60% glycerol fresh weekly.

Prepare the enrichment beads (macro-scale: 4 ePCR reactions)

1. Vortex the enrichment beads, then immediately transfer 825 μ L of the enrichment beads to each of two 2.0-mL LoBind Tubes.
2. Centrifuge the beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
3. Resuspend the enrichment beads in 500 μ L of 1 \times Bind & Wash Buffer per tube.
4. Combine the contents of the two tubes into a single tube, resulting in one 2.0-mL tube containing enrichment beads in 1 mL of 1 \times Bind & Wash Buffer.
5. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
6. Resuspend the enrichment beads in 500 μ L of 1 \times Bind & Wash Buffer.
7. Add 5 μ L of 1 mM Enrichment Oligo, then vortex and pulse-spin the enrichment beads.
8. Rotate the tube at room temperature for 30 minutes.
9. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
10. Resuspend the enrichment beads in 1 mL of 1 \times TEX Buffer.
11. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
12. Repeat steps 10 and 11.
13. Resuspend the enrichment beads in 500 μ L of 1 \times Low Salt Binding Buffer.

STOPPING POINT. Store the prepared enrichment beads at 4 °C in 1 \times Low Salt Binding Buffer, or proceed to “[Prepare the templated beads for enrichment \(macro-scale: 4 ePCR reactions\)](#)”. Prepared enrichment beads should be used within one week of preparation.

Prepare the templated beads for enrichment (macro-scale: 4 ePCR reactions)

1. Place a 2.0-mL LoBind Tube in a magnetic rack.
2. Transfer the suspension of templated beads from the first ePCR reaction to the tube in the magnetic rack.
3. Rinse the bottom of the first tube of templated beads with 100 μ L of 1 \times TEX Buffer, then transfer the rinse to the tube in the magnetic rack.
4. Wait for at least 1 minute. After the solution clears, remove and discard the supernatant.
5. Transfer the suspension of templated beads from the next ePCR reaction to the tube in the magnetic rack.
6. Rinse the tube with 100 μ L of 1 \times TEX Buffer, then transfer the rinse to the tube in the magnetic rack.
7. Wait for at least 1 minute. After the solution clears, remove and discard the supernatant.
8. Repeat steps 5 to 7 until all templated beads are in the LoBind Tube in the magnetic rack.
9. Resuspend the templated beads in 450 μ L of prepared Denaturing Buffer solution, then let the mixture stand for 1 minute.
10. Place the tube in a magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
11. Repeat steps 9 and 10 twice.
12. Resuspend the beads in 1.0 mL of 1 \times TEX Buffer.
13. Place the tube in a magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
14. Repeat steps 12 and 13 twice.
15. Resuspend the beads in 500 μ L of 1 \times TEX Buffer.
16. Sonicate the beads with the Covalent Declump 1 program on the Covaris[™] S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)).

Enrich the templated beads (macro-scale: 4 ePCR reactions)

1. Transfer all (500 μ L) of the enrichment bead suspension to the 2.0-mL tube with the templated beads, vortex to mix, then pulse-spin the tube.
2. Sonicate the enrichment-templated bead mixture using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 3” on page 127](#)), then pulse-spin the beads.
3. Incubate the bead mixture at 61 °C for 15 minutes. During the incubation, vortex and pulse-spin the bead mixture every 5 minutes including at the end of the incubation.
4. Immediately cool the beads on ice for 2 minutes.
5. Add 10 mL of *freshly prepared* 60% glycerol to a new 50-mL conical polypropylene tube.
6. Use a 1-mL pipettor tip to pipet the bead mixture up and down to mix, then load the entire volume of bead mixture *carefully* on top of the 60% glycerol solution. Do *not* vortex the tube.
7. Centrifuge the tube for 10 minutes at $3400 \times g$ (minimum $2284 \times g$). The centrifuge brake should be off and the temperature should be set to room temperature.
8. Add 10 mL of 1× TEX Buffer to a new 50-mL conical polypropylene tube.
9. Transfer the top layer of beads into the tube with 1× TEX Buffer. Aspirate as little glycerol as possible to collect all of the beads at the top layer without touching the un-templated beads at the bottom of the tube. When you dispense the top layer of beads into the 1× TEX Buffer, dispense the beads into the bottom of the tube. Aspirate a small amount of 1× TEX buffer to clean the pipette tip.
10. Top off the tube with additional 1× TEX Buffer to the 25-mL mark, then vortex the tube.
11. Centrifuge the tube for 10 minutes at $3400 \times g$ (minimum $2284 \times g$).



Note: Verify that the beads are pelleted in case excess glycerol carried over to the 1× TEX Buffer creates a matrix that impedes pelleting of beads.

12. Proceed according to the [Table 16](#):

Table 16 Steps for pelleted or unpelleted beads

If the beads are...	Then...
Pelleted	Remove and discard the supernatant, then proceed to “Isolate the P2-enriched beads (macro-scale: 4 ePCR reactions)” on page 80 .

Table 16 Steps for pelleted or unpelleted beads

If the beads are...	Then...
Not pelleted	Perform steps 13 to 15.

13. Carefully remove as much supernatant as possible without pipetting up the beads.
14. Top off the tube with additional 1× TEX Buffer to the 25-mL mark, then vortex the tube.
15. Repeat steps 11 to 12.

**Isolate the
P2-enriched beads
(macro-scale: 4
ePCR reactions)**

1. Resuspend the beads in 900 µL of prepared Denaturing Buffer solution, then transfer the beads into a new 2.0-mL LoBind Tube. Let the beads stand for 1 minute.
2. Rinse the 50-mL tube with 300 µL of prepared Denaturing Buffer solution, then transfer the rinse to the same 2.0-mL LoBind Tube.
3. Place the tube in a magnetic rack for at least 1 minute until the supernatant is pure white or clear, then remove and discard the supernatant.

❗ **IMPORTANT!** *Never* magnet the P2-enriched beads before adding prepared Denaturing Buffer solution to the beads. If you do, the templated beads linked to the enrichment beads are lost when the supernatant is removed.
4. Resuspend the beads with 1 mL of prepared Denaturing Buffer solution, then let the beads stand for 1 minute.
5. Repeat steps 3 and 4 until the supernatant is clear (that is, until all white enrichment beads have been removed).
6. Place the tube in a magnetic rack for at least 1 minute until the supernatant is clear. Remove and discard the supernatant.
7. Resuspend the beads in 1 mL of 1× TEX Buffer.
8. Repeat steps 6 and 7 *twice*.
9. Sonicate the enrichment-templated bead mixture using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)), then pulse-spin the beads.
10. Place the tube in a magnetic rack for at least 1 minute until the supernatant is clear. Remove and discard the supernatant.
11. Resuspend the beads in 1 mL of 1× TEX Buffer.
12. Place the tube in a magnetic rack for at least 1 minute until the supernatant is clear. Remove and discard the supernatant.

13. If the supernatant appears cloudy due to residual enrichment beads, repeat steps 11 and 12 until the supernatant is clear.
14. Resuspend the beads in 1 mL of 1× TEX Buffer.

STOPPING POINT. Store the prepared enrichment beads at 4 °C in 1× TEX Buffer, or proceed to [“Modify the 3’ ends \(macro-scale: 4 ePCR reactions\)” on page 82.](#)

Modify the 3' ends (macro-scale: 4 ePCR reactions)

Extend the 3' ends with Terminal Transferase and Bead Linker (macro-scale: 4 ePCR reactions)

1. If the P2-enriched beads have been stored overnight or longer, sonicate the beads using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 3” on page 127](#)). Pulse-spin the beads.
2. Prepare the appropriate volume of 1× Terminal Transferase Reaction Buffer (1.5 mL per 4 ePCR reactions; see [Table 17](#)):

Table 17 Four ePCR reactions: prepare 1× Terminal Transferase Reaction Buffer

Component	Volume per reaction (μL)
10× Terminal Transferase Buffer	165
10× Cobalt Chloride	165
Nuclease-free Water	1170
Total	1500



Note: The 1× Terminal Transferase Reaction Buffer should be clear. If the solution becomes colored, discard it and then prepare a fresh buffer using a new lot of material.

3. Add 2 μL of 50 mM Bead Linker to 98 μL of 1× Low TE Buffer to prepare a 1 mM Bead Linker solution.
4. Place the tube of P2-enriched beads in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
5. Resuspend the beads in 300 μL of 1× Terminal Transferase Reaction Buffer, then transfer the beads to a new 2.0-mL LoBind Tube.
6. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
7. Resuspend the beads in 300 μL of 1× Terminal Transferase Reaction Buffer.
8. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
9. Resuspend the beads in 712 μL of 1× Terminal Transferase Reaction Buffer.
10. Add 80 μL of 1 mM Bead Linker solution to the tube.
11. Sonicate the beads using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 3” on page 127](#)). Pulse-spin the beads.

12. Add 8.0 μL of Terminal Transferase (20 U/ μL) to the tube, vortex, then pulse-spin the beads.
13. Seal the tube with Parafilm, then place the tube on a rotator and rotate for 2 hours at 37 °C.
14. Pulse-spin the tube.
15. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
16. Resuspend the beads in 400 μL of 1 \times TEX Buffer.
17. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
18. Resuspend the beads in 400 μL of 1 \times TEX Buffer.

STOPPING POINT. Store the templated beads at 4 °C in 1 \times TEX Buffer, or proceed to “Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer (macro-scale: 4 ePCR reactions)” or “Quantitate the beads using a hemocytometer” on page 115.

Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer (macro-scale: 4 ePCR reactions)

1. If necessary, generate a standard curve (see “Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer” on page 112).
2. Sonicate the beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see “Covalent Declump 1” on page 127), then pulse-spin the beads.
3. Use the SOLiD™ Bead Concentration Chart (Applied Biosystems PN 4415131) to estimate the bead concentration of the beads (see Figure 37 for a picture of the chart; for best results use the official chart).

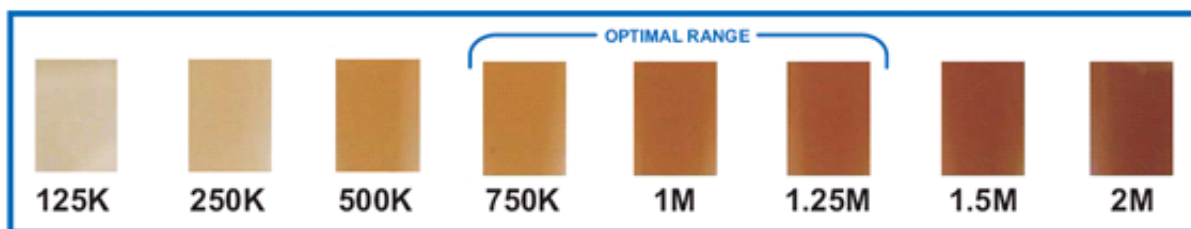


Figure 37 The SOLiD™ Bead Concentration Chart (facsimile). For best results, use the SOLiD™ Bead Concentration Chart (PN 4415131), supplied separately.

4. Adjust the volume of beads so that the color of the bead solution matches a color in the optimal range (750,000 beads/ μL to 1.25 million beads/ μL ; see Figure 38 on page 84 for the workflow).

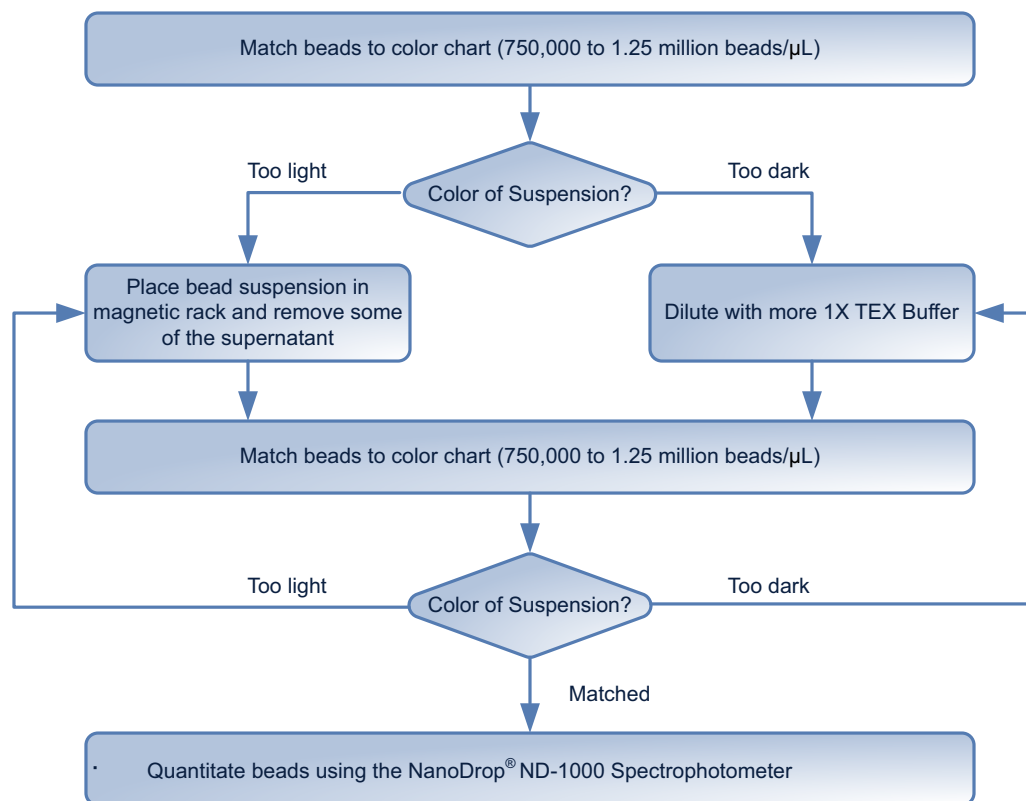


Figure 38 The SOLiD™ Bead Concentration Chart workflow.

5. When the bead concentration is within accurate range, quantitate the beads using the NanoDrop® ND-1000 Spectrophotometer. Take three readings, then average them. Calculate the bead concentration using the appropriate standard curve (for more details, see [“Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer”](#) on page 112).

STOPPING POINT. Store the templated beads at 4 °C in 1X TEX Buffer, or proceed to bead deposition and sequencing [refer to the *Applied Biosystems SOLiD™ 4 System Instrument Operation Guide* (PN 4448379)].

7

Prepare Macro-Scale Templated Beads (8 ePCR Reactions)

- Prepare the templated beads (macro-scale: 8 ePCR reactions). 86
- Prepare 4 full-scale emulsion PCR (ePCR) reactions. 86
- Perform the emulsion breaks and bead wash (macro-scale: 8 ePCR reactions). . 86
- Enrich the templated beads (macro-scale: 8 ePCR reactions) 86
- Modify the 3' ends (macro-scale: 8 ePCR reactions) 91

Prepare the templated beads (macro-scale: 8 ePCR reactions)

These instructions describe how to generate 1.2 billion to 2.4 billion templated beads using the macro-scale (8 ePCR reactions) templated bead preparation method.

Note: Do not begin these instructions unless you have completed the instructions in [Chapter 5, Macro-Scale Templated Beads \(4 and 8 ePCR Reactions\)](#). [Chapter 5](#) also describes the workflow overview.

To instead prepare macro-scale (4 ePCR reactions), use [Chapter 6, “Prepare Macro-Scale Templated Beads \(4 ePCR Reactions\)”](#) on page 75.

Prepare 4 full-scale emulsion PCR (ePCR) reactions

Prepare 8 emulsion PCR reactions as described in “[Prepare the full-scale emulsion PCR \(ePCR\) reaction](#)” in [Chapter 4, Prepare Full-Scale Templated Beads](#). Eight ePCR reactions provide an adequate bead yield for one full slide for sequencing. Store each 96-well plate at 4 °C or proceed to “[Perform the emulsion breaks and bead wash \(macro-scale: 8 ePCR reactions\)](#)”.


Perform the emulsion breaks and bead wash (macro-scale: 8 ePCR reactions)

Perform the emulsion break and bead wash procedure on each of the 8 emulsion PCR reactions, as described in “[Perform the emulsion break and bead wash \(full-scale\)](#)” in [Chapter 4, Prepare Full-Scale Templated Beads](#). Store each tube of beads at 4 °C or proceed to “[Enrich the templated beads \(macro-scale: 8 ePCR reactions\)](#)”.

Enrich the templated beads (macro-scale: 8 ePCR reactions)

Prepare the Denaturing Buffer solution


1. For each enrichment (8 plates to be combined), transfer 8.1 mL of Denaturing Buffer to a 15-mL conical tube.
2. Add 900 µL of Denaturant to the 8.1 mL of Denaturing Buffer, then cap the tube and vortex.

 **IMPORTANT!** Create a new batch of the prepared Denaturing Buffer solution each week.

Prepare 60% glycerol

1. With a 10-mL syringe, add 8 mL of Nuclease-free Water to a 50-mL conical tube.

2. With a 3-mL syringe, add 12 mL of glycerol to the Nuclease-free Water by dispensing 3 mL of glycerol four times with the syringe. Fill and dispense the glycerol slowly to ensure that the total volume of glycerol is dispensed.
3. Cap the tube, then vortex to mix the solution well.

 **IMPORTANT!** Prepare a new solution of 60% glycerol each week.

**Prepare the
enrichment beads
(macro-scale: 8
ePCR reactions)**

1. Vortex the enrichment beads, then immediately transfer 825 μ L of the enrichment beads to each of four 2.0-mL LoBind Tubes.
2. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
3. Resuspend the enrichment beads in 500 μ L of 1 \times Bind & Wash Buffer per tube.
4. Combine the contents of two tubes into a single tube, resulting in two 2.0-mL tubes containing enrichment beads, each tube with 1 mL of 1 \times Bind & Wash Buffer.
5. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
6. Resuspend the enrichment beads in 500 μ L of 1 \times Bind & Wash Buffer per tube.
7. Add 5 μ L of 1 mM Enrichment Oligo *per tube*, then vortex and pulse-spin the tube of enrichment beads.
8. Rotate the tubes at room temperature for 30 minutes.
9. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
10. Resuspend the enrichment beads in 1 mL of 1 \times TEX Buffer per tube.
11. Centrifuge the enrichment beads for 2 minutes at $21,000 \times g$ (minimum $14,000 \times g$), then remove and discard the supernatant.
12. Repeat steps 10 and 11.
13. Resuspend the enrichment beads in each of the two tubes with 500 μ L of 1 \times Low Salt Binding Buffer per tube.

STOPPING POINT. Store the prepared enrichment beads at 4 °C in 1 \times Low Salt Binding Buffer, or proceed to “[Prepare the templated beads for enrichment \(macro-scale: 4 ePCR reactions\)](#)” on page 78. Prepared enrichment beads should be used within one week of preparation.

Prepare the templated beads for enrichment (macro-scale: 8 ePCR reactions)

1. Place a 2.0-mL LoBind Tube in a magnetic rack.
2. Transfer the suspension of templated beads from the first ePCR reaction to the tube in the magnetic rack.
3. Rinse the bottom of the first tube of templated beads with 100 μ L of 1 \times TEX Buffer, then transfer the rinse to the tube in the magnetic rack.
4. Wait for at least 1 minute. After the solution clears, remove and discard the supernatant.
5. Transfer the suspension of templated beads from the next ePCR reaction to the tube in the magnetic rack.
6. Rinse the tube with 100 μ L of 1 \times TEX Buffer, then transfer the rinse to the tube in the magnetic rack.
7. Wait for at least 1 minute. After the solution clears, remove and discard the supernatant.
8. Repeat steps 5 to 7 until all templated beads from four tubes are in the LoBind Tube in the magnetic rack.
9. Repeat steps 1 to 8 for the remaining four tubes of templated beads.
10. Resuspend the templated beads in each tube with 450 μ L of prepared Denaturing Buffer solution, then let stand for 1 minute.
11. Place the tubes in a magnetic rack for at least 1 minute. After the solution clears, remove then discard the supernatant.
12. Repeat steps 10 and 11 *twice*.
13. Resuspend the beads in each tube with 1.0 mL of 1 \times TEX Buffer.
14. Place the tubes in a magnetic rack for at least 1 minute. After the solution clears, remove and discard the supernatant.
15. Repeat steps 13 and 14 *twice*.
16. Resuspend the beads in each tube with 500 μ L of 1 \times TEX Buffer.
17. Sonicate the beads with the Covalent Declump 1 program on the Covaris[™] S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)).

**Enrich the
 templated beads
 (macro-scale: 8
 ePCR reactions)**

1. Transfer one tube of enrichment beads (500 μ L) into one 2.0-mL tube of templated beads. Vortex to mix, then pulse spin the tube.
2. Immediately sonicate the enrichment-templated bead mixture using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)), then pulse-spin the beads.
3. Repeat steps 1 and 2 for the second tube of enrichment and templated beads. You will have 2 tubes of enrichment and templated bead mixtures, each containing 1 mL of total volume.
4. Incubate each tube of the bead mixture at 61 °C for 15 minutes, then vortex and pulse-spin the bead mixture every 5 minutes including at the end of the incubation.
5. Immediately cool the beads on ice for 2 minutes.
6. Add 10 mL of *freshly prepared* 60% glycerol to each of two new 50-mL conical polypropylene tubes.
7. Use a 1-mL pipettor tip to pipet the bead mixture from one of the tubes up and down to mix, then load the entire volume of bead mixture *carefully* on top of one of the 60% glycerol solutions. *Do not* vortex the tube.
8. Centrifuge the tubes for 10 minutes at $3400 \times g$ (minimum $2284 \times g$). The centrifuge brake should be off and the temperature should be set to room temperature.
9. Repeat step 8 for the second tube of enrichment and templated bead mixture and the second tube of 60% glycerol solution.
10. Add 20 mL of 1× TEX Buffer to a new 50-mL conical polypropylene tube.
11. Transfer the top layer of beads from each glycerol cushion into the tube with 1× TEX Buffer. Aspirate as little glycerol as possible to collect all of the beads at the top layer without touching the un-templated beads at the bottom of the tube. When you dispense the top layer of beads into the 1× TEX Buffer, dispense the beads into the bottom of the tube. Aspirate a small amount of 1× TEX buffer to clean the pipette tip.
12. Top off the tube with additional 1× TEX Buffer to the 35-mL mark, then vortex the tube.

13. Centrifuge the tube for 10 minutes at $3400 \times g$ (minimum $2284 \times g$).



Note: Verify that the beads are pelleted in case excess glycerol carried over to the 1× TEX Buffer creates a matrix that impedes pelleting of beads.

14. Proceed according to [Table 18](#):

Table 18 Steps for pelleted or unpelleted beads

If the beads are...	Then perform steps...
Pelleted	Remove and discard the supernatant, then proceed to “Isolate the P2-enriched beads (macro-scale: 8 ePCR reactions)” on page 90.
Not pelleted	Perform steps 13 to 15.

15. Carefully remove as much supernatant as possible without pipetting up the beads.
16. Top off the tube with additional 1× TEX Buffer to the 25-mL mark, then vortex the tube.
17. Repeat steps 13 and 14.

Isolate the P2-enriched beads (macro-scale: 8 ePCR reactions)

1. Resuspend the beads in 900 μL of prepared Denaturing Buffer solution, then transfer the beads into a new 2.0-mL LoBind Tube. Let the beads stand for 1 minute.
2. Rinse the 50-mL tube with 300 μL of prepared Denaturing Buffer solution, then transfer the rinse to the same 2.0-mL LoBind Tube.
3. Place the tube in a magnetic rack for at least 1 minute until the supernatant is pure white or clear, then remove and discard the supernatant.

⚠ **IMPORTANT!** *Never* magnet the P2-enriched beads before adding prepared Denaturing Buffer solution to the beads. If you do, the templated beads linked to the enrichment beads are lost when the supernatant is removed.
4. Resuspend the beads with 1 mL of prepared Denaturing Buffer solution, then let the beads stand for 1 minute.
5. Repeat steps 3 and 4 until the supernatant is clear (all white enrichment beads have been removed).

6. Place the tube in a magnetic rack for at least 1 minute until the supernatant is clear. Remove and discard the supernatant.
7. Resuspend the beads in 1 mL of 1× TEX Buffer.
8. Repeat steps 6 and 7 *twice*.
9. Sonicate the enrichment-templated bead mixture using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)), then pulse-spin the beads.
10. Place the tube in a magnetic rack for at least 1 minute until the supernatant is clear. Remove and discard the supernatant.
11. Resuspend the beads in 1 mL of 1× TEX Buffer.
12. Place the tube in a magnetic rack for at least 1 minute until the supernatant is clear. Remove and discard the supernatant.
13. If the supernatant appears cloudy due to residual enrichment beads, repeat steps 11 and 12 until the supernatant is clear.
14. Resuspend the beads in 1 mL of 1× TEX Buffer.

STOPPING POINT. Store the prepared enrichment beads at 4 °C in 1× TEX Buffer, or proceed to [“Modify the 3' ends \(macro-scale: 8 ePCR reactions\)”](#).

Modify the 3' ends (macro-scale: 8 ePCR reactions)

Extend the 3' ends with Terminal Transferase and Bead Linker

1. If the P2-enriched beads have been stored overnight or longer, sonicate the beads using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)). Pulse-spin the beads.
2. Prepare the appropriate volume of 1× Terminal Transferase Reaction Buffer (2.4 mL per 8 ePCR reactions; see [Table 19](#)).

Table 19 Eight ePCR reactions: prepare 1× Terminal Transferase Reaction Buffer

Component	Volume per reaction (μL)
10× Terminal Transferase Buffer	264
10× Cobalt Chloride	264
Nuclease-free Water	1872
Total	2400



Note: The 1× Terminal Transferase Reaction Buffer should be clear. If the solution becomes colored, discard then prepare fresh buffer using a new lot of material.

3. Add 4 μL of 50 mM Bead Linker to 196 μL of 1× Low TE Buffer to prepare a 1 mM Bead Linker solution.
4. Place the tube of P2-enriched beads in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
5. Resuspend the beads in 300 μL of 1× Terminal Transferase Reaction Buffer, then transfer the beads to a 2.0-mL LoBind Tube.
6. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
7. Resuspend the beads in 300 μL of 1× Terminal Transferase Reaction Buffer.
8. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
9. Resuspend the beads in 1424 μL of 1× Terminal Transferase Reaction Buffer.
10. Add 160 μL of 1 mM Bead Linker solution to the tube.
11. Transfer 792 μL of bead solution to a new 2.0-mL LoBind Tube.
12. Sonicate the beads using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 1” on page 127](#)). Pulse-spin the beads.
13. Add 8 μL of Terminal Transferase (20 U/μL) to each tube, vortex, then pulse-spin the beads.
14. Seal the tubes with Parafilm, then place the tubes on a rotator and rotate for 2 hours at 37 °C.
15. Pulse-spin the tubes, then pool the beads in one LoBind Tube.

16. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
17. Resuspend the beads in 400 μ L of 1 \times TEX Buffer.
18. Place the tube in a magnetic rack for at least 1 minute. After the supernatant clears, remove and discard the supernatant.
19. Resuspend the beads in 400 μ L of 1 \times TEX Buffer.

STOPPING POINT. Store the templated beads at 4 $^{\circ}$ C in 1 \times TEX Buffer, or proceed to “Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer (macro-scale: 8 ePCR reactions)” or “Quantitate the beads using a hemocytometer” on page 115.

Quantitate the beads with the SOLiD™ Bead Concentration Chart and the NanoDrop® ND-1000 Spectrophotometer (macro-scale: 8 ePCR reactions)

1. If necessary, generate a standard curve (see “Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer” on page 112).
2. Sonicate the beads using the Covalent Declump 1 program on the Covaris™ S2 System (for program conditions, see “Covalent Declump 1” on page 127), then pulse-spin the beads.
3. Use the SOLiD™ Bead Concentration Chart (Applied Biosystems PN 4415131) to estimate the bead concentration of the beads (see Figure 39 for a picture of the chart; for best results use the official chart).

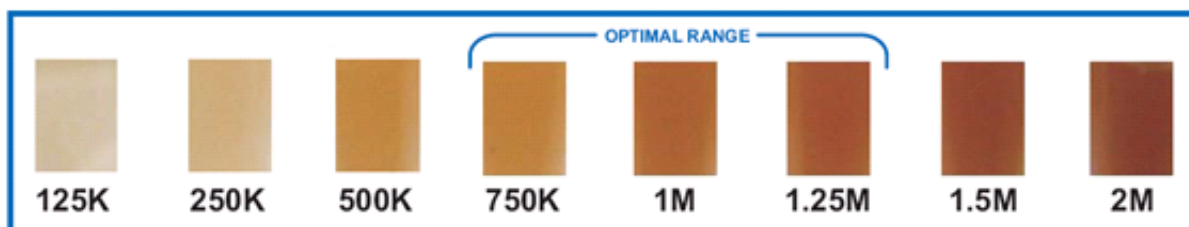


Figure 39 The SOLiD™ Bead Concentration Chart (facsimile). For best results, use the SOLiD™ Bead Concentration Chart (PN 4415131), supplied separately.

4. Adjust the volume of beads so that the color of the bead solution matches a color in the optimal range (750,000 beads/ μ L to 1.25 million beads/ μ L; see Figure 40 for the workflow).

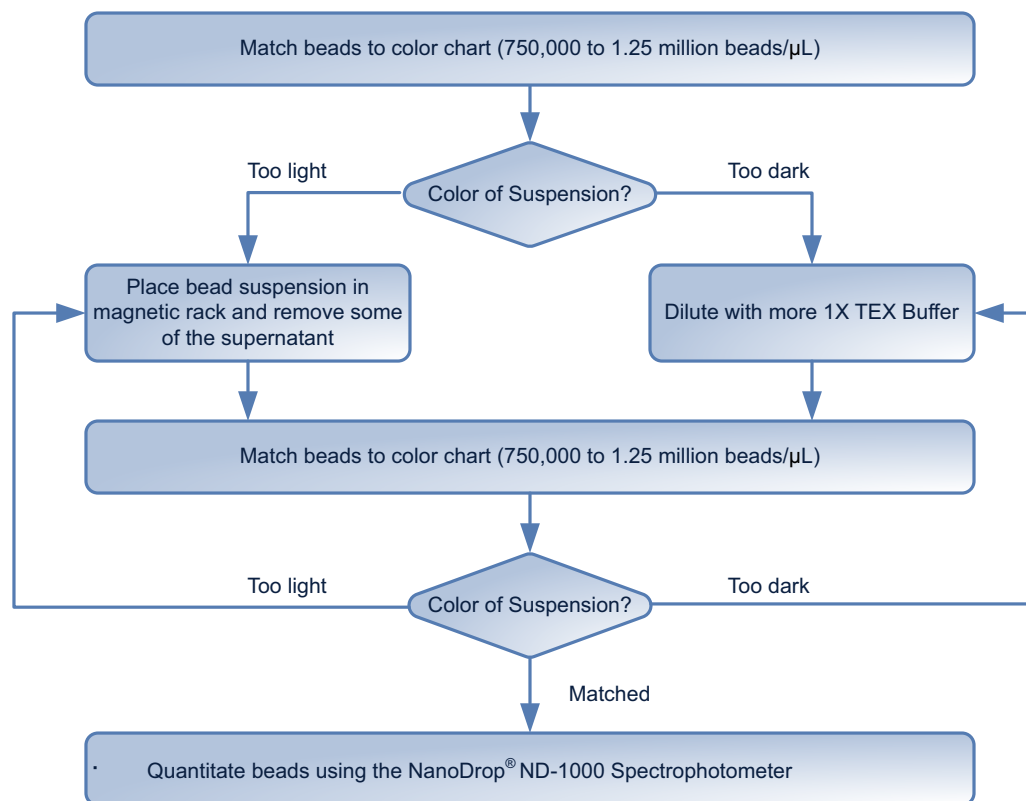


Figure 40 The SOLiD™ Bead Concentration Chart workflow.

5. When the bead concentration is within accurate range, quantitate the beads using the NanoDrop® ND-1000 Spectrophotometer. Take three readings, then average them. Calculate the bead concentration using the appropriate standard curve (for more details, see [“Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer”](#) on page 112).

STOPPING POINT. Store the templated beads at 4 °C in 1X TEX Buffer, or proceed to bead deposition and sequencing [refer to the *Applied Biosystems SOLiD™ 4 System Instrument Operation Guide* (PN 4448379)].



Required Materials

- Prepare templated beads (mini-scale) 96
- Prepare templated beads (full-scale and macro-scale) 100

Prepare templated beads (mini-scale)

Table 20 Required Applied Biosystems reagent kits

Item (part number) [‡]	Components	Kit component(s) used in...
SOLiD™ ePCR Kit V2, 20 Mini-Reactions (4407756)	Magnesium Chloride Emulsion Oil Emulsion Stabilizer 1 Emulsion Stabilizer 2 Bead Block Solution 10X PCR Buffer dNTP Mix AmpliTaq Gold DNA Polymerase, UP ePCR primer 1 ePCR primer 2 SOLiD™ P1 DNA Beads	Emulsion PCR
SOLiD™ Buffer Kit, 20 Mini-Reactions (4407759)	1X Bead Wash Buffer 2-Butanol [§]	Emulsion break and bead wash
	1X Bind & Wash Buffer 1X Low Salt Binding Buffer	Enrichment
	1X Low TE Buffer	Emulsion PCR, 3'-end modification
	1X TEX Buffer	Emulsion PCR, emulsion break and bead wash, enrichment, 3'-end modification
SOLiD™ Bead Enrichment Kit, 20 Mini-Reactions # (4407757)	Glycerol Denaturing Buffer Denaturant Enrichment Oligo Enrichment Beads	Enrichment
SOLiD™ Pre Deposition Kit (4445808)	10X Terminal Transferase Buffer 10X Cobalt Chloride Terminal Transferase Bead Linker	3'-end modification

[‡] Applied Biosystems has validated this protocol using this specific material. Substitution may adversely affect system performance.

[§] The tube is labeled as “butanol” in the kit.

[#] If more Denaturing Buffer and Denaturant are needed, order the SOLiD™ Bead Enrichment Kit Box 1 of 3 (4392177), which contains Denaturing Buffer, Denaturant, and Glycerol.

Table 21 Required equipment

Item [‡]	Source
ULTRA-TURRAX® Tube Drive from IKA®§ (115 V for U.S. customers) (230 V for international customers) The system includes: SOLiD™™ ePCR Tubes and Caps, 10-pack	Applied Biosystems 4400335 (115 V and 230 V)
96-well GeneAmp® PCR System 9700 (thermal cycler)	<ul style="list-style-type: none"> Applied Biosystems N8050200 (Base) Applied Biosystems 4314443 (Block)[‡]
Covaris™ S2 System (110 V for U.S. customers) (220 V for international customers) The system includes: <ul style="list-style-type: none"> Covaris™ S2 sonicator Latitude™ laptop from Dell® Inc. MultiTemp III Thermostatic Circulator Covaris-2 series Machine Holder for (one) 1.5-mL microcentrifuge tube Covaris-2 series Machine Holder for (one) 0.65-mL microcentrifuge tube Covaris-2 series Machine Holder for (one) 13 mm × 65 mm tube Covaris-2 Series Machine Holder for (one) microTUBE Covaris microTUBE Prep Station Covaris Water Tank Label Kit Covaris microTUBES (1 pack of 25) For the materials summary for the Covaris™ S2 System, refer to the <i>Applied Biosystems SOLiD™ Site Preparation Guide</i> .	<ul style="list-style-type: none"> Applied Biosystems 4387833 (110 V) Applied Biosystems 4392718 (220 V) or Covaris
6-Tube Magnetic Stand	Applied Biosystems AM10055
Microcentrifuge 5417R, refrigerated, without rotor	<ul style="list-style-type: none"> Eppendorf[#] 022621807 (120 V/60 Hz) Eppendorf[‡] 022621840 (230 V/50 Hz)
FA-45-24-11, fixed-angle rotor, 24 × 1.5/2 mL, including aluminum lid, aerosol-tight	Eppendorf [#] 022636006
Repeater® Xstream	Eppendorf 022460811

Table 21 Required equipment (continued)

Item [‡]	Source
Repeater [®] Plus Pipette	Eppendorf 022260201
NanoDrop [®] ND-1000 Spectrophotometer (computer required)	Thermo Scientific ND-1000
Labquake Rotisserie Rotator, Barnstead/Thermolyne	Thermo Scientific 400110
Fume hood	Major Laboratory Supplier (MLS)
Tabletop Centrifuge	MLS
Vortexer	MLS
Picofuge	MLS
Incubator (37 °C)	MLS
Incubator (61 °C)	MLS
Pipettors, 2 µL	MLS
Pipettors, 20 µL	MLS
Pipettors, 200 µL	MLS
Pipettors, 1000 µL	MLS

‡ Applied Biosystems has validated this protocol using this specific material. Substitution may adversely affect system performance.

§ Applied Biosystems ships one ULTRA-TURRAX[®] Tube Drive from IKA[®] per instrument.

In some cases equivalent equipment may be substituted. Validation of the equipment for library preparation is required.

Table 22 Required consumables

Item [‡]	Source
SOLiD [™] ePCR Tubes and Caps, 10 pack (15-mL tubes)	Applied Biosystems 4400401
SOLiD [™] Emulsion Collection Tray Kit	Applied Biosystems 4415129
MicroAmp [®] Optical 96-Well Reaction Plates	Applied Biosystems N8010560
Clear Adhesive Film: MicroAmp [®] Optical Adhesive Film, or Clear Seal Diamond Heat Sealing Film	MicroAmp [®] Optical Adhesive Film: Applied Biosystems 4360954 Clear Seal Diamond Heat Sealing Film: Thermo Scientific, AB-0812
Nuclease-free Water (1 L)	Applied Biosystems AM9932
50-mL high-clarity polypropylene conical centrifuge tube, 9400 RCF rating, sterile	Becton-Dickinson 352070

Table 22 Required consumables (continued)

Item [‡]	Source
1-mL BD™ slip-tip disposable tuberculin syringe	Becton-Dickinson 309602
5-mL Combitips Plus	Eppendorf 022496107
10-mL Combitips Plus	Eppendorf 022496123
0.5-mL LoBind Tubes	Eppendorf 022431005
1.5-mL LoBind Tubes	Eppendorf 022431021
2.0-mL LoBind Tubes	Eppendorf 022431048
Polypropylene wide-mouth jars (0.5 oz., 15 mL, 38-mm cap)	Nalgene 2118-9050
Ethylene glycol	American Bioanalytical AB00455-01000
CF-1 Calibration Fluid Kit	Thermo Scientific CF-1
PR-1 Conditioning Kit [§]	Thermo Scientific PR-1
10-mL serological pipettes	Major Laboratory Supplier (MLS)
15-mL conical polypropylene tubes	MLS
3-mL syringes	MLS
10-mL syringes	MLS
Tape	MLS
Razor blades	MLS
Filtered pipettor tips	MLS
Ice	MLS

[‡] Applied Biosystems has validated this protocol using this specific material. Substitution may adversely affect system performance.

[§] The NanoDrop® Conditioning Kit is useful for reconditioning the sample measurement pedestals to a hydrophobic state if they become unconditioned (refer to the NanoDrop™ user's manual for more information). The PR-1 kit consists of a container of specially formulated polishing compound and a supply of convenient applicators.

Prepare templated beads (full-scale and macro-scale)

Table 23 Required Applied Biosystems reagent kits

Item (part number) [‡]	Components	Kit component(s) used in...
SOLiD™ ePCR Kit V2 (4400834)	Magnesium Chloride Emulsion Oil Emulsion Stabilizer 1 Emulsion Stabilizer 2 Bead Block Solution 10× PCR Buffer dNTP Mix AmpliTaq Gold DNA Polymerase, UP ePCR primer 1 ePCR primer 2 SOLiD™ P1 DNA Beads	Emulsion PCR
SOLiD™ Buffer Kit (4387918)	1X Bead Wash Buffer 2-Butanol [§]	Emulsion break and bead wash
	1X Bind & Wash Buffer 1X Low Salt Binding Buffer	Enrichment
	1X Low TE Buffer	Emulsion PCR, 3'-end modification
	1X TEX Buffer	Emulsion PCR, emulsion break and bead wash, enrichment, 3'-end modification
SOLiD™™ XD Bead Enrichment Kit (4453663)	Glycerol Denaturing Buffer Denaturant Enrichment Oligo Enrichment Beads	Enrichment
SOLiD™™ Pre Deposition Kit (4445808)	10X Terminal Transferase Buffer 10X Cobalt Chloride Terminal Transferase Bead Linker	3'-end modification

[‡] Applied Biosystems has validated this protocol using this specific material. Substitution may adversely affect system performance.

[§] The tube is labeled as “butanol” in the kit.

Table 24 Required equipment

Item [‡]	Source
ULTRA-TURRAX [®] Tube Drive from IKA ^{®§} (115 V for U.S. customers) (230 V for international customers) The system includes: SOLiD [™] ePCR Tubes and Caps, 10-pack	Applied Biosystems 4400335 (115 V and 230 V)
96-well GeneAmp [®] PCR System 9700 (thermal cycler)	<ul style="list-style-type: none"> Applied Biosystems N8050200 (Base) Applied Biosystems 4314443 (Block)[‡]
Covaris [™] S2 System (110 V for U.S. customers) (220 V for international customers) The system includes: <ul style="list-style-type: none"> Covaris[™] S2 sonicator Latitude[™] laptop from Dell[®] Inc. MultiTemp III Thermostatic Circulator Covaris-2 series Machine Holder for (one) 1.5-mL microcentrifuge tube Covaris-2 series Machine Holder for (one) 0.65-mL microcentrifuge tube Covaris-2 series Machine Holder for (one) 13 mm × 65 mm tube Covaris-2 Series Machine Holder for (one) microTUBE Covaris microTUBE Prep Station Covaris Water Tank Label Kit Covaris microTUBEs (1 pack of 25) For the materials summary for the Covaris [™] S2 System, refer to the <i>Applied Biosystems SOLiD[™] Site Preparation Guide</i> .	<ul style="list-style-type: none"> Applied Biosystems 4387833 (110 V) Applied Biosystems 4392718 (220 V) or Covaris Inc.
6-Tube Magnetic Stand	Applied Biosystems AM10055
Microcentrifuge 5417R, refrigerated, without rotor	<ul style="list-style-type: none"> Eppendorf[#] 022621807 (120 V/60 Hz) Eppendorf[‡] 022621840 (230 V/50 Hz)
FA-45-24-11, fixed-angle rotor, 24 × 1.5/2 mL, including aluminum lid, aerosol-tight	Eppendorf [#] 022636006
Repeater [®] Xstream	Eppendorf 022460811

Table 24 Required equipment (continued)

Item [‡]	Source
Repeater [®] Plus Pipette	Eppendorf 022260201
NanoDrop [®] ND-1000 Spectrophotometer (computer required)	Thermo Scientific ND-1000
Labquake Rotisserie Rotator, Barnstead/Thermolyne	Thermo Scientific 400110
Fume hood	Major Laboratory Supplier (MLS)
Tabletop Centrifuge	MLS
Vortexer	MLS
Picofuge	MLS
Incubator (37 °C)	MLS
Incubator (61 °C)	MLS
Pipettors, 2 µL	MLS
Pipettors, 20 µL	MLS
Pipettors, 200 µL	MLS
Pipettors, 1000 µL	MLS

[‡] Applied Biosystems has validated this protocol using this specific material. Substitution may adversely affect system performance.

[§] Applied Biosystems ships one ULTRA-TURRAX[®] Tube Drive from IKA[®] per instrument.

[#] In some cases equivalent equipment may be substituted. Validation of the equipment for library preparation is required.

Table 25 Required consumables

Item [‡]	Source
SOLiD [™] ePCR Tubes and Caps, 10 pack (15-mL tubes)	Applied Biosystems 4400401
SOLiD [™] Emulsion Collection Tray Kit	Applied Biosystems 4415129
MicroAmp [®] Optical 96-Well Reaction Plates	Applied Biosystems N8010560
Clear Adhesive Film: MicroAmp [®] Optical Adhesive Film, or Clear Seal Diamond Heat Sealing Film	MicroAmp [®] Optical Adhesive Film: Applied Biosystems 4360954 Clear Seal Diamond Heat Sealing Film: Thermo Scientific, AB-0812
Nuclease-free Water (1 L)	Applied Biosystems AM9932
50-mL high-clarity polypropylene conical centrifuge tube, 9400 RCF rating, sterile	Becton-Dickinson 352070
1-mL BD [™] slip-tip disposable tuberculin syringe	Becton-Dickinson 309602

Table 25 Required consumables (continued)

Item [‡]	Source
5-mL Combitips Plus	Eppendorf 022496107
10-mL Combitips Plus	Eppendorf 022496123
0.5-mL LoBind Tubes	Eppendorf 022431005
1.5-mL LoBind Tubes	Eppendorf 022431021
2.0-mL LoBind Tubes	Eppendorf 022431048
Polypropylene wide-mouth jars (0.5 oz., 15 mL, 38-mm cap)	Nalgene 2118-9050
Ethylene glycol	American Bioanalytical AB00455-01000
CF-1 Calibration Fluid Kit	Thermo Scientific CF-1
PR-1 Conditioning Kit [§]	Thermo Scientific PR-1
10-mL serological pipettes	Major Laboratory Supplier (MLS)
15-mL conical polypropylene tubes	MLS
3-mL syringes	MLS
10-mL syringes	MLS
Tape	MLS
Razor blades	MLS
Filtered pipettor tips	MLS
Parafilm	MLS
Ice	MLS

[‡] Applied Biosystems has validated this protocol using this specific material. Substitution may adversely affect system performance.

[§] The NanoDrop[®] Conditioning Kit is useful for reconditioning the sample measurement pedestals to a hydrophobic state if they become unconditioned (refer to the Nanodrop[™] user's manual for more information). The PR-1 kit consists of a container of specially formulated polishing compound and a supply of convenient applicators.



Appendix A Required Materials

Prepare templated beads (full-scale and macro-scale)



Supplemental Procedures

- Program the Eppendorf Repeater® Xstream Pipettor 106
- Quantitate the beads using the NanoDrop® ND-1000 Spectrophotometer. 107
- Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer 112
- Quantitate the beads using a hemocytometer 115

Program the Eppendorf Repeater® Xstream Pipettor

The Eppendorf Repeater® Xstream pipettor has been preset to use with IKA®-based emulsions and the 10-mL Combitip Plus. Follow the procedure below only if you need to reprogram the pipettor.

Materials and equipment required

Table 26 Required equipment

Item	Source
Repeater® Xstream	Eppendorf 022460811

Table 27 Required consumables

Item	Source
10-mL Combitips Plus	Eppendorf 022496123

Procedure

1. Attach a 10-mL Combitip Plus on the Eppendorf Repeater Xstream pipettor.
2. Set the top dial to pipette mode: **Pip**.
3. Push the left blue **select** button. The screen displays “Set volume.”
4. Toggle the right blue +/- button to set the pipettor fill volume to **5.6 mL** (or other appropriate volume as specified in the procedure).
5. Push the left blue **select** button. The screen displays “up (▲) speed.”
6. Toggle the right blue +/- button to set histogram to **scale 5** (five colored bars: mid-range).
7. Push the left blue **select** button. The screen displays “down (▼) speed.”
8. Toggle the right blue +/- button to set histogram to **scale 1** (one colored bar: slowest).
9. Push the left blue **select** button to finish programming.
10. Push the *round lower center* blue button to save/store program.
11. Use the programmed Eppendorf Repeater Xstream pipettor with IKA®-based emulsions.

Quantitate the beads using the NanoDrop® ND-1000 Spectrophotometer

Materials and equipment required

Table 28 Required equipment

Item	Source
Covaris™ S2 System (110 V for U.S. customers) (220 V for international customers) The system includes: <ul style="list-style-type: none"> • Covaris™ S2 sonicator • Latitude™ laptop from Dell® Inc. • MultiTemp III Thermostatic Circulator • Covaris-2 series Machine Holder for (one) 1.5-mL microcentrifuge tube • Covaris-2 series Machine Holder for (one) 0.65-mL microcentrifuge tube • Covaris-2 series Machine Holder for (one) 13 mm × 65 mm tube • Covaris-2 Series Machine Holder for (one) microTUBE • Covaris microTUBE Prep Station • Covaris Water Tank Label Kit • Covaris microTUBEs (1 pack of 25) For the materials summary for the Covaris™ S2 System, refer to the <i>Applied Biosystems SOLiD™ Site Preparation Guide</i> .	Applied Biosystems PN 4387833 (110 V) PN 4392718 (220 V) or Covaris Inc.
NanoDrop® ND-1000 Spectrophotometer (computer required)	Thermo Scientific ND-1000
Pipettors	Major Laboratory Supplier (MLS)

Table 29 Required consumables

Item	Source
SOLiD™ Buffer Kit – 1X TEX Buffer	Applied Biosystems PN 4389776‡
Nuclease-free Water (1 L)	Applied Biosystems PN AM9932
CF-1 Calibration Fluid Kit§	Thermo Scientific CF-1
PR-1 Conditioning Kit	Thermo Scientific PR-1

Table 29 Required consumables

Item	Source
0.5-mL LoBind Tubes	Eppendorf 022431005
Filtered pipettor tips	Major Laboratory Supplier (MLS)

‡ The part number for the complete SOLiD™™ Buffer Kit is 4387918.

§ The NanoDrop® Conditioning Kit is useful for reconditioning the sample measurement pedestals to a hydrophobic state if they become unconditioned. (Refer to the NanoDrop user's manual for more information.) The PR-1 kit consists of a container of specially formulated polishing compound and a supply of convenient applicators.

Procedure

1. Ensure that the NanoDrop ND-1000 Spectrophotometer is properly calibrated. Use the CF-1 Calibration Fluid Kit if necessary.
2. Open the NanoDrop ND-1000 Spectrophotometer software. A dialog box displays (see [Figure 41](#)).

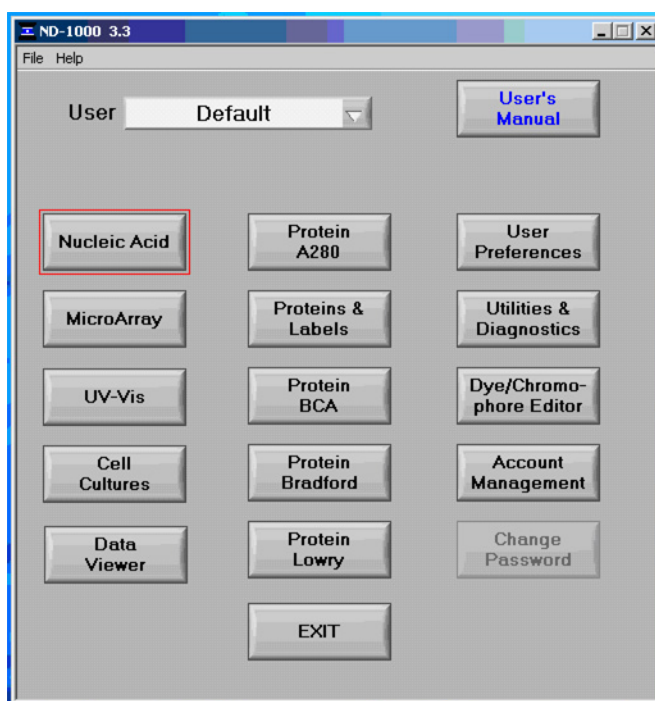


Figure 41 NanoDrop® ND-1000 Spectrophotometer software dialog box.

3. Select the **Cell Cultures** button.
4. Lift the sampling arm and load 2 µL of Nuclease-free Water onto the lower measurement pedestal and lower the sampling arm (see [Figure 42 on page 109](#)).

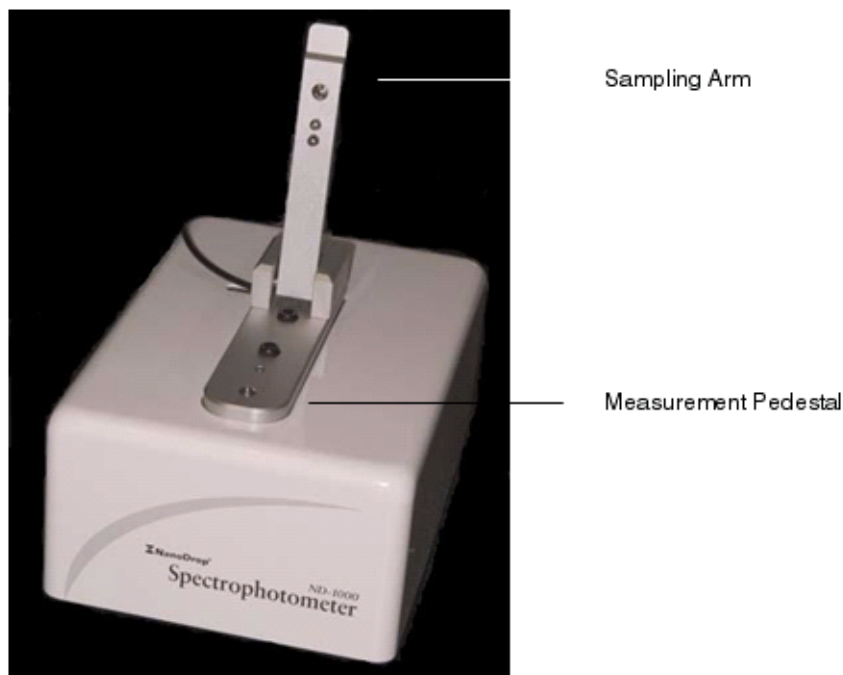


Figure 42 Components of the NanoDrop® ND-1000 Spectrophotometer.

5. In the dialog box, click **OK**, then allow the instrument to initialize.
6. Lift the sampling arm and use a Kimwipe® to remove water from the measurement pedestal and the sampling arm.
7. Load 2 μ L of the same buffer that was used to resuspend the beads onto the sampling pedestal, then lower the sampling arm.
8. Click **Blank**, then allow the instrument to take a measurement (see [Figure 43 on page 110](#)).

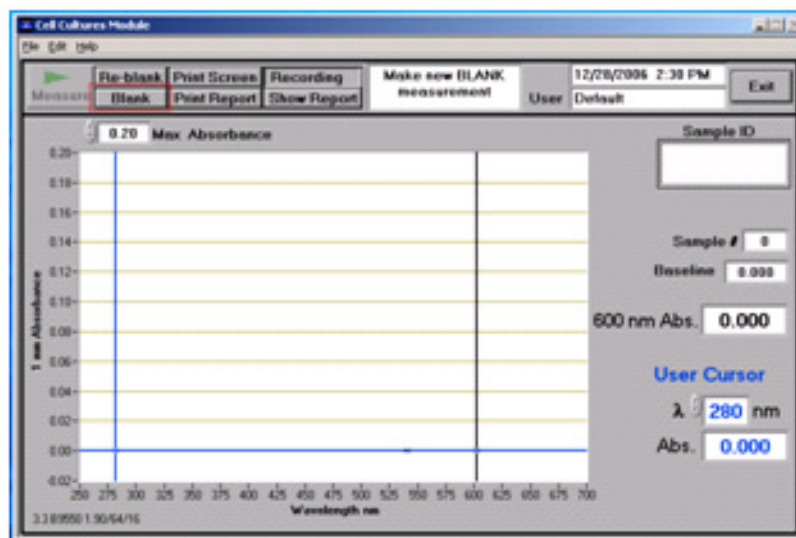


Figure 43 NanoDrop® ND-1000 Spectrophotometer software measurement dialog box.

9. Lift the sampling arm and wipe away the buffer from the sampling arm and measurement pedestal with a Kimwipe. The instrument is now ready to take readings.
10. Sonicate the beads using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 3” on page 127](#)), then pulse-spin the beads. Proceed immediately to the next step.
11. If necessary, make a dilution of beads in 1× TEX Buffer.
12. Lift the sampling arm and load 2 µL of beads onto the lower measurement pedestal and lower the sampling arm.
13. Enter the sample name in the Sample ID field and click **Measure**. The A600 readings should be between 0.2 and 1 absorbance unit. Depending on the absorbance, perform one of these steps:
 - If the absorbance reading is >1 abs, dilute beads until the absorbance reading is within the correct range.
 - If the absorbance reading is <0.2 abs, place the tube of beads in the magnetic rack and resuspend them in half the volume of buffer. Be sure to sonicate the beads again according to step 10.
14. Record the absorbance for each sample.
15. Use a Kimwipe® to clean the sample from the sampling arm and the measurement pedestal.
16. Repeat steps 12 to 15 two more times for a total of three readings.
17. Repeat steps 9 to 16 for any remaining samples.

18. (Optional) Save the data as a text document:
 - a. Click **Show Report** to open the Data Viewer.
 - b. Select **Reports ▶ Save Report As**.
 - c. Click the **Export Report Table Only** button to save the file in the desired location (see [Figure 44](#)).

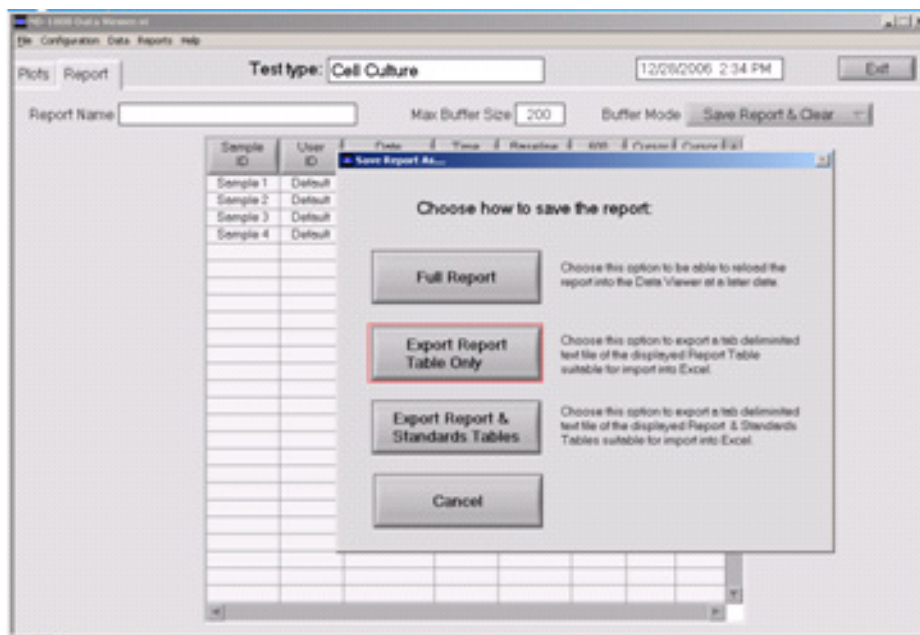


Figure 44 Save Report Software Dialog Box on the NanoDrop® ND-1000 Spectrophotometer.

19. Average the three A600 readings for each sample and calculate the bead concentrations using the appropriate standard curve (see [“Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer”](#) on page 112).

Generate a standard curve for calculating bead concentration using the NanoDrop® ND-1000 Spectrophotometer

Materials and equipment required

Table 30 Required equipment

Item	Source
The Covaris™ S2 System (110 V for U.S. customers) (220 V for international customers) The system includes: <ul style="list-style-type: none"> • Covaris™ S2 sonicator • Latitude™ laptop from Dell® Inc. • MultiTemp III Thermostatic Circulator • Covaris-2 series Machine Holder for (one) 1.5-mL microcentrifuge tube • Covaris-2 series Machine Holder for (one) 0.65-mL microcentrifuge tube • Covaris-2 series Machine Holder for (one) 13 mm × 65 mm tube • Covaris-2 Series Machine Holder for (one) microTUBE • Covaris microTUBE Prep Station • Covaris Water Tank Label Kit • Covaris microTUBEs (1 pack of 25) For system materials summary, refer to “Covaris™ S2 System Materials Summary,” in the <i>Applied Biosystems SOLiD™ Site Preparation Guide</i> .	Applied Biosystems PN 4387833 (110 V) PN 4392718 (220 V) or Covaris Inc.
NanoDrop® ND-1000 Spectrophotometer (computer required)	Thermo Scientific ND-1000
Hemocytometer	Major Laboratory Supplier (MLS)
Clicker counter	MLS
Pipettors	MLS

Table 31 Required consumables

Item	Source
SOLiD™ Buffer Kit – 1X TEX Buffer	Applied Biosystems PN 4389776‡
SOLiD™ ePCR Kit – P1 DNA Beads	Applied Biosystems PN 4392175§
0.5-mL LoBind Tubes	Eppendorf 022431005
Filtered pipettor tips	MLS

‡ The part number for the complete SOLiD™™ Buffer Kit is 4387918.

§ The part number for the complete SOLiD™™ ePCR Kit V2 is 4400834.

- Procedure**
1. Sonicate either P1 DNA Beads or surplus templated beads using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, [“Covalent Declump 3” on page 127](#)), then pulse-spin the beads.
 2. Dilute the beads to a concentration of between 10,000 and 100,000 beads/μL.
 3. Place the glass coverslip on the hemocytometer.
 4. Pipet 10 μL of diluted beads into the groove of the hemocytometer. Allow the beads to settle for 5 minutes.
 5. Count an average of 4 squares of the 25 squares that form the larger center square. Use a clicker counter and count the beads only within the triple lines of the square.
 6. Calculate the concentration of beads using the following formula:
Bead concentration = (average beads in square) × 250 × (dilution factor)
 - Example**
Bead concentration = (240 beads) × 250 × 100 = 6.0×10^6 beads/μL
 7. Rinse, then dry the hemocytometer.
 8. According to the hemocytometer counts, dilute ePCR beads in 1× TEX to make 10 μL of the following concentrations: 200 K, 400 K, 600 K, 800 K, 1 M, and 1.2 M beads/μL, where K = 10^3 and M = 10^6 .
 9. Take readings on the NanoDrop® ND-1000 Spectrophotometer for each bead concentration (see [“Quantitate the beads using the NanoDrop® ND-1000 Spectrophotometer” on page 107](#)). The lowest absorbance reading should be < 0.2 and the largest absorbance readings should be > 1. If the above dilution series does not meet these criteria, create additional dilutions).
 10. Using analysis software such as Microsoft® Office Excel®, average the NanoDrop readings for each concentration and graphically plot absorbance versus bead concentration. A linear trend line gives the equation of the standard curve, $y = mx + b$ (see Figure 4), where:
y: Absorbance at 600 nm
m: Slope of the line
x: Bead concentration (beads/μL)
b: y-intercept (determined by extrapolating standard curve) (see [Figure 45 on page 114](#))

Examples

SOLiD™ P1 DNA Beads: $(A_{600}) = (8 \times 10^{-7}) \times (\text{Concentration}) + 0.08$

Enrichment Beads: $(A_{600}) = (4 \times 10^{-6}) \times (\text{Concentration}) + 0.04$

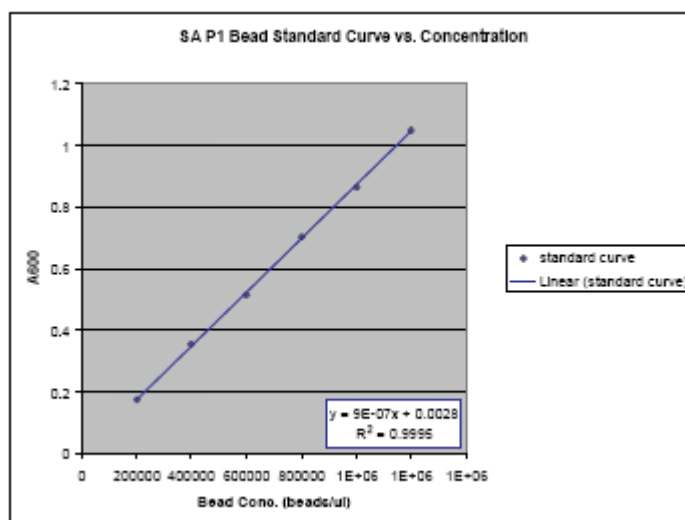


Figure 45 Standard curve generated from NanoDrop® readings of a titration of beads.

11. For added accuracy, repeat steps 1 to 9 with new dilutions and average the resulting curves.
12. Create an Excel analysis worksheet to convert a NanoDrop reading to concentration. The formula is:

$$\text{Concentration} = \frac{(\text{A600 Measurement}) - (y\text{-intercept})}{\text{slope}}$$

Quantitate the beads using a hemocytometer

Materials and equipment required

Table 32 Required equipment

Item	Source
Covaris™ S2 System (110 V for U.S. customers) (220 V for international customers) The system includes: <ul style="list-style-type: none"> • Covaris™ S2 sonicator • Latitude™ laptop from Dell® Inc. • MultiTemp III Thermostatic Circulator • Covaris-2 series Machine Holder for (one) 1.5-mL microcentrifuge tube • Covaris-2 series Machine Holder for (one) 0.65-mL microcentrifuge tube • Covaris-2 series Machine Holder for (one) 13 mm × 65 mm tube • Covaris-2 Series Machine Holder for (one) microTUBE • Covaris microTUBE Prep Station • Covaris Water Tank Label Kit • Covaris microTUBEs (1 pack of 25) For the materials summary for the Covaris™ S2 System, refer to the <i>Applied Biosystems SOLiD™ Site Preparation Guide</i> .	Applied Biosystems PN 4387833 (110 V) PN 4392718 (220 V) or Covaris Inc.
Hemocytometer	Major Laboratory Supplier (MLS)
Microscope	MLS
Clicker counter	MLS
Pipettors	MLS [‡]

[‡] For the Safety Data Sheet (SDS) of any chemical not distributed by Applied Biosystems, contact the chemical manufacturer. Before handling any chemicals, refer to the SDS provided by the manufacturer, and observe all relevant precautions.

Table 33 Required consumables

Item	Source
SOLiD™ Buffer Kit – 1X TEX Buffer	Applied Biosystems PN 4389776 [‡]
0.5-mL LoBind Tubes	Eppendorf 022431005
Filtered pipettor tips	MLS

[‡] The part number for the complete SOLiD™ Buffer Kit is 4387918.

- Procedure**
1. Sonicate the beads using the Covalent Declump 3 program on the Covaris™ S2 System (for program conditions, see [“Covalent Declump 3” on page 127](#)), then pulse-spin the beads.
 2. Make a dilution of beads in 1× TEX Buffer (1:100 dilution recommended for post-emulsion break quantitation and 1:10 dilution recommended for post-3' -end-modification quantitation).
 3. Place the glass coverslip on the hemocytometer.
 4. Pipet 10 µL of diluted beads into the groove of the hemocytometer. Allow the beads to settle for 5 minutes.
 5. Count beads in 4 squares of the 25 squares that form the larger center square. Use a clicker counter and count the beads only within the triple lines of the square. Determine the average of your 4 counts.
 6. Calculate the concentration of beads using the following formula:
Bead concentration = (average beads in square) × 250 × (dilution factor)
Example
Bead concentration = (240 beads) × 250 × 100 = 6.0×10^6 beads/µL
 7. Rinse, then dry the hemocytometer.

C

Library Concentration Conversion

Fragment Library:

Assuming average length of 215 bp

$$X \text{ pg/}\mu\text{L DNA} = \frac{500 \text{ pmol}}{1 \text{ L}} \times \frac{660 \text{ pg}}{1 \text{ pmol}} \times \frac{1 \text{ L}}{10^6 \mu\text{L}} \times 215 = 71 \text{ pg/}\mu\text{L DNA}$$

2 × 50 bp Mate-Paired Library:

Assuming average length of 300 bp

$$X \text{ pg/}\mu\text{L DNA} = \frac{500 \text{ pmol}}{1 \text{ L}} \times \frac{660 \text{ pg}}{1 \text{ pmol}} \times \frac{1 \text{ L}}{10^6 \mu\text{L}} \times 300 = 99 \text{ pg/}\mu\text{L DNA}$$

2 × 25 bp Mate-Paired Library:

Assuming average length of 155 bp

$$X \text{ pg/}\mu\text{L DNA} = \frac{500 \text{ pmol}}{1 \text{ L}} \times \frac{660 \text{ pg}}{1 \text{ pmol}} \times \frac{1 \text{ L}}{10^6 \mu\text{L}} \times 155 = 51 \text{ pg/}\mu\text{L DNA}$$

Barcoded Fragment Library:

Assuming average length of 255 bp

$$X \text{ pg/}\mu\text{L DNA} = \frac{500 \text{ pmol}}{1 \text{ L}} \times \frac{660 \text{ pg}}{1 \text{ pmol}} \times \frac{1 \text{ L}}{10^6 \mu\text{L}} \times 255 = 84 \text{ pg/}\mu\text{L DNA}$$

Whole Transcriptome Library:

Assuming average length of 230 bp

$$X \text{ pg/}\mu\text{L DNA} = \frac{500 \text{ pmol}}{1 \text{ L}} \times \frac{660 \text{ pg}}{1 \text{ pmol}} \times \frac{1 \text{ L}}{10^6 \mu\text{L}} \times 175 = 76 \text{ pg/}\mu\text{L DNA}$$

Small RNA Library:

Assuming average length of 123 bp

$$X \text{ pg/}\mu\text{L DNA} = \frac{500 \text{ pmol}}{1 \text{ L}} \times \frac{660 \text{ pg}}{1 \text{ pmol}} \times \frac{1 \text{ L}}{10^6 \mu\text{L}} \times 130 = 41 \text{ pg/}\mu\text{L DNA}$$

Barcoded SAGE Library:

Assuming average length of 130 bp

$$X \text{ pg/}\mu\text{L DNA} = \frac{500 \text{ pmol}}{1 \text{ L}} \times \frac{660 \text{ pg}}{1 \text{ pmol}} \times \frac{1 \text{ L}}{10^6 \mu\text{L}} \times 130 = 43 \text{ pg/}\mu\text{L DNA}$$





Checklists and Workflow Tracking Forms

- Workflow checklists: prepare templated beads. 120
- Workflow tracking: prepare templated beads (mini-scale or full-scale). 122
- Workflow tracking: prepare templated beads (macro-scale) 123

Workflow checklists: prepare templated beads

	Equipment	Reagents	Preparation steps
Emulsion PCR (ePCR)	<input type="checkbox"/> Covaris™ S2 System <input type="checkbox"/> ULTRA-TURRAX® Tube Drive from IKA® <input type="checkbox"/> SOLiD™ ePCR Tube <input type="checkbox"/> Thermal cycler <input type="checkbox"/> Xstream Pipettor <input type="checkbox"/> Repeater Plus Pipette <input type="checkbox"/> Magnetic rack <input type="checkbox"/> Vortexer <input type="checkbox"/> Picofuge <input type="checkbox"/> 3-mL syringe <input type="checkbox"/> 1-mL syringe <input type="checkbox"/> 10-mL syringe <input type="checkbox"/> 5-mL Combitip Plus <input type="checkbox"/> 10-mL Combitip Plus <input type="checkbox"/> 1.5-mL LoBind tubes <input type="checkbox"/> 50-mL conical polypropylene tubes <input type="checkbox"/> Pipettors <input type="checkbox"/> Filtered pipettor tips <input type="checkbox"/> 96-well PCR plates <input type="checkbox"/> Wide-mouthed jars <input type="checkbox"/> Clear adhesive film <input type="checkbox"/> Razor blades <input type="checkbox"/> Ice	<input type="checkbox"/> Library template <input type="checkbox"/> SOLiD™ P1 DNA Beads <input type="checkbox"/> Emulsion Stabilizer 1 <input type="checkbox"/> Emulsion Stabilizer 2 <input type="checkbox"/> Emulsion Oil <input type="checkbox"/> ePCR Primer 1 <input type="checkbox"/> ePCR Primer 2 <input type="checkbox"/> 10× PCR Buffer <input type="checkbox"/> dNTP Mix <input type="checkbox"/> Magnesium chloride <input type="checkbox"/> 1× Low TE Buffer <input type="checkbox"/> 1× TEX Buffer <input type="checkbox"/> Bead Block Solution <input type="checkbox"/> Nuclease-free water	<input type="checkbox"/> Turn on Covaris™ S2 System (including chiller and degasser) <input type="checkbox"/> Thaw library template, ePCR Primer 1, ePCR Primer 2, dNTP Mix, 10× PCR Buffer
Emulsion break and bead wash	<input type="checkbox"/> Covaris™ S2 System <input type="checkbox"/> NanoDrop® ND-1000 <input type="checkbox"/> Fume hood <input type="checkbox"/> Tabletop centrifuge <input type="checkbox"/> Microcentrifuge <input type="checkbox"/> Magnetic rack <input type="checkbox"/> Vortexer <input type="checkbox"/> Picofuge <input type="checkbox"/> SOLiD™ Emulsion Collection Tray <input type="checkbox"/> Tape <input type="checkbox"/> 50-mL conical polypropylene tubes <input type="checkbox"/> 1.5-mL LoBind tubes <input type="checkbox"/> Pipettors <input type="checkbox"/> Filtered pipettor tips <input type="checkbox"/> Paper towels	<input type="checkbox"/> 2-butanol <input type="checkbox"/> 1× Bead Wash Buffer <input type="checkbox"/> 1× TEX Buffer	<input type="checkbox"/> Turn on Covaris™ S2 System (including chiller and degasser)

	Equipment	Reagents	Preparation steps
Templated bead enrichment	<input type="checkbox"/> Covaris™ S2 System <input type="checkbox"/> Incubator (61 °C) <input type="checkbox"/> Microcentrifuge <input type="checkbox"/> Rotator <input type="checkbox"/> Magnetic rack <input type="checkbox"/> Vortexer <input type="checkbox"/> PicoFuge <input type="checkbox"/> 0.5-mL Lo-Bind tube <input type="checkbox"/> 1.5-mL Lo-Bind tube <input type="checkbox"/> 2.0-mL Lo-Bind tube <input type="checkbox"/> 3-mL syringe <input type="checkbox"/> 10-mL syringe <input type="checkbox"/> 15-mL conical polypropylene tubes <input type="checkbox"/> Pipettors <input type="checkbox"/> Filtered pipettor tips <input type="checkbox"/> Ice	<input type="checkbox"/> Enrichment Beads <input type="checkbox"/> Enrichment Oligo <input type="checkbox"/> Denaturing Buffer <input type="checkbox"/> Denaturant <input type="checkbox"/> Glycerol <input type="checkbox"/> 1× Bind & Wash Buffer <input type="checkbox"/> 1× TEX Buffer <input type="checkbox"/> 1× Low Salt Binding Buffer <input type="checkbox"/> Nuclease-free water	<input type="checkbox"/> Turn on Covaris™ S2 System (including chiller and degasser) <input type="checkbox"/> Turn on 61 °C incubator <input type="checkbox"/> Thaw Enrichment Oligo
3' -End modification	<input type="checkbox"/> Covaris™ S2 System <input type="checkbox"/> NanoDrop® ND-1000 <input type="checkbox"/> Incubator (37 °C) <input type="checkbox"/> Rotator <input type="checkbox"/> Magnetic rack <input type="checkbox"/> Vortexer <input type="checkbox"/> PicoFuge <input type="checkbox"/> 1.5-mL LoBind tubes <input type="checkbox"/> Pipettors <input type="checkbox"/> Filtered pipettor tips	<input type="checkbox"/> 10× Terminal Transferase Buffer <input type="checkbox"/> 10× Cobalt chloride <input type="checkbox"/> Terminal Transferase <input type="checkbox"/> Bead Linker <input type="checkbox"/> 1× Low TE Buffer <input type="checkbox"/> 1× TEX Buffer <input type="checkbox"/> Nuclease-free water	<input type="checkbox"/> Turn on Covaris™ S2 System (including chiller and degasser) <input type="checkbox"/> Turn on 37 °C incubator <input type="checkbox"/> Thaw 10× Terminal Transferase Buffer and Bead Linker

Workflow tracking: prepare templated beads (mini-scale or full-scale)

Sample:			
Quantitation			
After emulsion break & bead wash		After 3' -end modification	
Lot numbers			
Emulsion PCR (ePCR)		Templated bead enrichment	
SOLiD™ ePCR Kit Box 1 of 3		SOLiD™ Bead Enrichment Kit Box 1 of 3	
SOLiD™ ePCR Kit Box 2 of 3		SOLiD™ Bead Enrichment Kit Box 2 of 3	
SOLiD™ ePCR Kit Box 3 of 3		SOLiD™ Bead Enrichment Kit Box 3 of 3	
Emulsion Stabilizer 1		Denaturing Buffer	
Emulsion Stabilizer 2		Denaturant	
Emulsion Oil		Glycerol	
SOLiD™ ePCR Tube		Enrichment Beads	
ePCR Primer 1		Enrichment Oligo	
ePCR Primer 2		1× Bind & Wash Buffer	
1× Low TE Buffer		1× TEX Buffer	
10× PCR Buffer		1× Low Salt Binding Buffer	
dNTP Mix		3' -End modification	
Magnesium chloride		SOLiD™ Pre Deposition Kit	
AmpliTaq Gold® DNA Polymerase, UP		10× Terminal Transferase Buffer	
SOLiD™ P1 DNA Beads		10× Cobalt chloride	
Bead Block Solution		Bead Linker	
1× TEX Buffer		Terminal Transferase	
Emulsion break and bead wash		1× TEX Buffer	
2-butanol			
SOLiD™ Emulsion Collection Tray			
1× Bead Wash Buffer			
1× TEX Buffer			

Sample:			
Quantitation			
After emulsion break & bead wash		After 3' -end modification	
Lot numbers			
Emulsion PCR (ePCR)		Templated bead enrichment	
SOLiD™ ePCR Kit Box 1 of 3		SOLiD™ Bead Enrichment Kit Box 1 of 3	
SOLiD™ ePCR Kit Box 2 of 3		SOLiD™ Bead Enrichment Kit Box 2 of 3	
SOLiD™ ePCR Kit Box 3 of 3		SOLiD™ Bead Enrichment Kit Box 3 of 3	
Emulsion Stabilizer 1		Denaturing Buffer	
Emulsion Stabilizer 2		Denaturant	
Emulsion Oil		Glycerol	
SOLiD™ ePCR Tube		Enrichment Beads	
ePCR Primer 1		Enrichment Oligo	
ePCR Primer 2		1× Bind & Wash Buffer	
1× Low TE Buffer		1× TEX Buffer	
10× PCR Buffer		1× Low Salt Binding Buffer	
dNTP Mix		3' -End modification	
Magnesium chloride		SOLiD™ Pre Deposition Kit	
AmpliTaq Gold® DNA Polymerase, UP		10× Terminal Transferase Buffer	
SOLiD™ P1 DNA Beads		10× Cobalt chloride	
Bead Block Solution		Bead Linker	
1× TEX Buffer		Terminal Transferase	
Emulsion break and bead wash		1× TEX Buffer	
2-butanol			
SOLiD™ Emulsion Collection Tray			
1× Bead Wash Buffer			
1× TEX Buffer			

Workflow tracking: prepare templated beads (macro-scale)

Sample:			
Quantitation			
After emulsion break & bead wash (Plate 1)		After 3' -end modification	
After emulsion break & bead wash (Plate 2)			
After emulsion break & bead wash (Plate 3)			
After emulsion break & bead wash (Plate 4)			
After emulsion break & bead wash (Plate 5)			
After emulsion break & bead wash (Plate 6)			
After emulsion break & bead wash (Plate 7)			
After emulsion break & bead wash (Plate 8)			
Lot numbers			
Emulsion PCR (ePCR)		Templated bead enrichment	
SOLiD™ ePCR Kit Box 1 of 3		SOLiD™ Bead Enrichment Kit Box 1 of 3	
SOLiD™ ePCR Kit Box 2 of 3		SOLiD™ Bead Enrichment Kit Box 2 of 3	
SOLiD™ ePCR Kit Box 3 of 3		SOLiD™ Bead Enrichment Kit Box 3 of 3	
Emulsion Stabilizer 1		Denaturing Buffer	
Emulsion Stabilizer 2		Denaturant	
Emulsion Oil		Glycerol	
SOLiD™ ePCR Tube		Enrichment Beads	
ePCR Primer 1		Enrichment Oligo	
ePCR Primer 2		1× Bind & Wash Buffer	
1× Low TE Buffer		1× TEX Buffer	
10× PCR Buffer		1× Low Salt Binding Buffer	
dNTP Mix		3' -End modification	
Magnesium chloride		SOLiD™ Pre Deposition Kit	
AmpliTaq Gold DNA Polymerase, UP		10× Terminal Transferase Buffer	
SOLiD™ P1 DNA Beads		10× Cobalt chloride	
Bead Block Solution		Bead Linker	
1× TEX Buffer		Terminal Transferase	
Emulsion break and bead wash		1× TEX Buffer	
2-butanol			
SOLiD™ Emulsion Collection Tray			
1× Bead Wash Buffer			
1× TEX Buffer			



Appendix D Checklists and Workflow Tracking Forms
Workflow tracking: prepare templated beads (macro-scale)



The Covaris™ S2 System

- Operation notes 126
 - Fill the tank 126
 - Degas the water 126
 - Set the chiller 126
 - Perform required maintenance of the Covaris™ S2 System 126
- Covaris™ S2 programs 127
 - Bead Block Declump 127
 - Covalent Declump 1 127
 - Covalent Declump 3 127

Operation notes



Note: For important instrument safety information, refer to the Covaris™ S2 System manual.

Fill the tank Fill the tank with fresh deionized water to the proper fill line (fill-line level “15” is recommended). The water should cover the visible part of the tube.

Degas the water Degas the water for 30 minutes. To maintain degassed water, keep the pump continuously on during operation and sample processing.

Set the chiller Set the chiller temperature to between 2 to 5 °C to ensure that the temperature reading in the water bath displays 5 °C. The circulated water chiller should be supplemented with 20% ethylene glycol.

Perform required maintenance of the Covaris™ S2 System The Covaris™ S2 System requires regular maintenance to work properly. Perform the tasks in the table below (see [Table 34](#)):

Table 34 Required maintenance of the Covaris™ S2 System

Required maintenance task	Frequency to perform task
Degas water for 30 minutes prior to use	Before every use
Change water	Daily
Clean with bleach	Every two weeks

Covaris™ S2 programs

Bead Block
Declump**Table 35** Bead Block Declump: 1 cycle Treatment 1 followed by 1 cycle Treatment 2

	Treatment 1	Treatment 2
Duty Cycle	1%	5%
Intensity	5	5
Cycles/Burst	50	100
Time	5 sec	60 sec
Target wattage power performance estimate (W) [‡]	2	10

[‡] Not programmedCovalent
Declump 1**Table 36** Covalent Declump 1: 1 cycle Treatment 1 followed by 1 cycle Treatment 2

	Treatment 1	Treatment 2
Duty Cycle	2%	5%
Intensity	6	9
Cycles/Burst	100	100
Time	5 sec	30 sec
Target wattage power performance estimate (W) [‡]	4	15

[‡] Not programmedCovalent
Declump 3**Table 37** Covalent Declump 3: 3 cycles Treatment 1 followed by 1 cycle Treatment 2

	Treatment 1	Treatment 2
Duty Cycle	2%	5%
Intensity	6	9
Cycles/Burst	100	100
Time	5 sec	30 sec
Target wattage power performance estimate (W) [‡]	4	15

[‡] Not programmed





Instrument Warranty Information

- Computer configuration 130
- Limited product warranty 130
- Warranty period effective date 131
- Warranty claims 131
- Warranty exceptions 131
- Warranty limitations 132
- Damages, claims, and returns 133

Computer configuration

Applied Biosystems supplies or recommends certain configurations of computer hardware, software, and peripherals for use with its instrumentation. Applied Biosystems reserves the right to decline support for or impose extra charges for supporting nonstandard computer configurations or components that have not been supplied or recommended by Applied Biosystems. Applied Biosystems also reserves the right to require that computer hardware and software be restored to the standard configuration prior to providing service or technical support. For systems that have built-in computers or processing units, installing unauthorized hardware or software may void the Warranty or Service Plan.

Limited product warranty

Applied Biosystems warrants that all standard components of the SOLiD™ 4 Analyzer, IKA® ULTRA-TURRAX® Tube Drive, the Covaris™ S2 System, APC UPS, and the recirculating chiller will be free of defects in materials and workmanship for a period of one (1) year from the date the warranty period begins. Applied Biosystems will repair or replace, at its discretion, all defective components during this warranty period. Applied Biosystems warrants the Genomic Solutions HydroShear® DNA Shearing Device will be free of defects in materials and workmanship for a period of one (1) year from the date the warranty period begins. Applied Biosystems will replace a defective HydroShear DNA Shearing Device during the warranty period. The following parts of the HydroShear DNA Shearing Device are use- replaceable and not covered by the warranty on the HydroShearDNA Shearing Device: shearing assembly, syringes, syringe adapters, syringe shields, and output tubing. Applied Biosystems reserves the right to use new, repaired, or refurbished instruments or components for warranty and post-warranty service agreement replacements. Repair or replacement of products or components that are under warranty does not extend the original warranty period.

Applied Biosystems warrants that all optional accessories supplied with its SOLiD™ 4 Analyzer, such as peripherals, printers, and special monitors, will be free of defects in materials and workmanship for a period of ninety (90) days from the date the warranty begins. Applied Biosystems will repair or replace, at its discretion, defective accessories during this warranty period. After this warranty period, Applied Biosystems will pass on to the buyer, to the extent that it is permitted to do so, the warranty of the original manufacturer for such accessories.

With the exception of consumable and maintenance items, replaceable products or components used on or in the instrument are themselves warranted to be free of defects in materials and workmanship for a period of ninety (90) days.

Applied Biosystems warrants that chemicals and other consumable products will be free of defects in materials and workmanship when received by the buyer, but not thereafter, unless otherwise specified in documentation accompanying the product.

Applied Biosystems warrants that for a period of ninety (90) days from the date the warranty period begins, the tapes, diskettes, or other media bearing the operating software of the product, if any, will be free of defects in materials and workmanship under normal use. If there is a defect in the media covered by the above warranty and the media is returned to Applied Biosystems within the ninety (90) day warranty period, Applied Biosystems will replace the defective media.

Unless indicated herein, Applied Biosystems makes no warranty whatsoever in regard to products or parts furnished by third parties, including but not limited to the non-APC- branded UPS or APC UPS, Covaris™ S2, Genomic Solutions HydroShear DNA Shearing Device, Recirculating Chiller, and IKA ULTRA-TURRAX purchased or obtained from a third party. Such products or parts will be subject to the warranties, if any, of their respective manufacturers to the extent they are 'transferable or otherwise available to Applied Biosystems' buyer.

Applied Biosystems at its sole discretion may refuse to provide buyer with support or service for buyer's use of Covaris™ S2 in a method not described in a SOLiD™ System protocol.

Applied Biosystems does not warrant that the operation of the instrument or its operating software will be uninterrupted or be error-free.

Warranty period effective date

Any applicable warranty period under these sections begins on the earlier of the date of installation or ninety (90) days from the date of shipment for hardware and software installed by Applied Biosystems personnel. For all hardware and software installed by the buyer or anyone other than Applied Biosystems, and for all other products, the applicable warranty period begins the date the product is delivered to the buyer.

Warranty claims

Warranty claims must be made within the applicable warranty period, or, for chemicals or other consumable products, within thirty (30) days after receipt by the buyer unless otherwise specified in the documentation accompanying the product.

Warranty exceptions

The above warranties do not apply to defects resulting from misuse, neglect, or accident, including without limitation: operation with incompatible solvents or samples in the system; operation outside of the environmental or use specifications or not in conformance with the instructions for the instrument system, software, or accessories; improper or inadequate maintenance by the user; installation of software or interfacing, or use in combination with software or products, not supplied or authorized by Applied Biosystems; modification or repair of the product not authorized by Applied

Biosystems; relocation or movement of the instrument by buyer or by any third party not acting on behalf of Applied Biosystems; or intrusive activity, including without limitation, computer viruses, hackers or other unauthorized interactions with instrument or software that detrimentally affects normal operations.

Parts in contact with any liquid are considered wetted and may be deemed user-replaceable and not be covered by the above warranties, including, but not limited to, seals, filters, gaskets, shearing assemblies, valves, syringes, syringe adapters, syringe shields, and output tubing.

Warranty limitations

THE FOREGOING PROVISIONS SET FORTH APPLIED BIOSYSTEMS' SOLE AND EXCLUSIVE REPRESENTATIONS, WARRANTIES, AND OBLIGATIONS WITH RESPECT TO THE PRODUCTS WARRANTED HEREIN, AND APPLIED BIOSYSTEMS MAKES NO OTHER WARRANTY OF ANY KIND WHATSOEVER, EXPRESSED OR IMPLIED, INCLUDING WITHOUT LIMITATION, WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE, WHETHER ARISING FROM A STATUTE OR OTHERWISE IN LAW OR FROM A COURSE OF DEALING OR USAGE OF TRADE, ALL OF WHICH ARE EXPRESSLY DISCLAIMED.

THE REMEDIES PROVIDED HEREIN ARE THE BUYER'S SOLE AND EXCLUSIVE REMEDIES. WITHOUT LIMITING THE GENERALITY OF THE FOREGOING, TO THE FULL EXTENT ALLOWED BY LAW, IN NO EVENT SHALL APPLIED BIOSYSTEMS BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE (INCLUDING WITHOUT LIMITATION, ANY TRADE PRACTICE, UNFAIR COMPETITION, OR OTHER STATUTE OF SIMILAR IMPORT) OR ON ANY OTHER BASIS, FOR DIRECT, INDIRECT, PUNITIVE, INCIDENTAL, MULTIPLE, CONSEQUENTIAL, OR SPECIAL DAMAGES SUSTAINED BY THE BUYER OR ANY OTHER PERSON OR ENTITY, WHETHER OR NOT FORESEEABLE AND WHETHER OR NOT APPLIED BIOSYSTEMS IS ADVISED OF THE POSSIBILITY OF SUCH DAMAGES, INCLUDING WITHOUT LIMITATION, DAMAGES ARISING FROM OR RELATED TO LOSS OF USE, LOSS OF DATA, FAILURE OR INTERRUPTION IN THE OPERATION OF ANY EQUIPMENT OR SOFTWARE, DELAY IN REPAIR OR REPLACEMENT, OR FOR LOSS OF REVENUE OR PROFITS, LOSS OF GOOD WILL, LOSS OF BUSINESS, OR OTHER FINANCIAL LOSS OR PERSONAL INJURY OR PROPERTY DAMAGE.

NO AGENT, EMPLOYEE, OR REPRESENTATIVE OF APPLIED BIOSYSTEMS HAS ANY AUTHORITY TO MODIFY THE TERMS OF THIS LIMITED WARRANTY STATEMENT OR TO BIND APPLIED BIOSYSTEMS TO ANY AFFIRMATION, REPRESENTATION, OR WARRANTY CONCERNING THE PRODUCT THAT IS NOT CONTAINED IN THIS LIMITED WARRANTY STATEMENT, AND ANY SUCH MODIFICATION, AFFIRMATION, REPRESENTATION, OR WARRANTY MADE BY ANY AGENT, EMPLOYEE, OR REPRESENTATIVE OF APPLIED BIOSYSTEMS WILL NOT BE BINDING ON APPLIED BIOSYSTEMS, UNLESS IN A WRITING SIGNED BY AN EXECUTIVE OFFICER OF APPLIED BIOSYSTEMS.

THIS WARRANTY IS LIMITED TO THE BUYER OF THE PRODUCT FROM APPLIED BIOSYSTEMS AND IS NOT TRANSFERABLE.

Some countries or jurisdictions limit the scope of or preclude limitations or exclusion of warranties, of liability, such as liability for gross negligence or willful misconduct, or of remedies or damages, as or to the extent set forth above. In such countries and jurisdictions, the limitation or exclusion of warranties, liability, remedies or damages set forth above shall apply to the fullest extent permitted by law, and shall not apply to the extent prohibited by law.

Damages, claims, and returns

- Damages** If shipping damage to the product is discovered, contact the shipping carrier and request inspection by a local agent. Secure a written report of the findings to support any claim. Do not return damaged goods to Applied Biosystems without first securing an inspection report and contacting Applied Biosystems Technical Support for a Return Authorization (RA) number.
- Claims** After a damage inspection report is received by Applied Biosystems, Applied Biosystems will process the claim unless other instructions are provided.
- Returns** Do not return any material without prior notification and authorization.
- If for any reason it becomes necessary to return material to Applied Biosystems, contact Applied Biosystems Technical Support or your nearest Applied Biosystems subsidiary or distributor for a return authorization (RA) number and forwarding address. Place the RA number in a prominent location on the outside of the shipping container, and return the material to the address designated by the Applied Biosystems representative.





Safety

■ Instrumentation safety	136
General instrument safety	136
Physical hazard safety	136
■ Chemical safety	137
General chemical safety	137
SDSs	138
Chemical waste safety	139
Biological hazard safety	141



Instrumentation safety

General instrument safety



Note: For important instrument safety information, refer to the *Applied Biosystems SOLiD™ 4 System Instrument Operation Guide* (PN 4448379) and the Covaris™ S2 System manual. For general safety information, see the “Preface” on page 7.

Operating the instrument

Ensure that anyone who operates the instrument has:

- Received instructions in both general safety practices for laboratories and specific safety practices for the instrument.
- Read and understood all applicable Safety Data Sheets (SDSs, formerly known as Material Safety Data Sheets). See “About SDSs” on page 138.

Cleaning or decontaminating the instrument



CAUTION! Using a cleaning or decontamination method other than that specified by the manufacturer may result in damage to the instrument.

Physical hazard safety

Solvents and pressurized fluids



WARNING! PHYSICAL INJURY HAZARD. Always wear eye protection when working with solvents or any pressurized fluids.

- Be aware that PEEK™ tubing is a polymeric material. Use caution when working with any polymer tubing that is under pressure.
Always wear eye protection when near pressurized polymer tubing.
- Extinguish all nearby flames if you use flammable solvents.
- Do not use PEEK tubing that has been severely stressed or kinked.
- Do not use PEEK tubing with tetrahydrofuran or nitric and sulfuric acids.
- Be aware that methylene chloride and dimethyl sulfoxide cause PEEK tubing to swell and greatly reduce the rupture pressure of the tubing.
- Be aware that high solvent flow rates (~40 mL/min) may cause a static charge to build up on the surface of the tubing. Electrical sparks may result.



Chemical safety

General chemical safety

Chemical hazard warning



WARNING! CHEMICAL HAZARD. Before handling any chemicals, refer to the Safety Data Sheet (SDS) provided by the manufacturer, and observe all relevant precautions.



WARNING! CHEMICAL HAZARD. All chemicals in the instrument, including liquid in the lines, are potentially hazardous. Always determine what chemicals have been used in the instrument before changing reagents or instrument components. Wear appropriate eyewear, protective clothing, and gloves when working on the instrument.



WARNING! CHEMICAL HAZARD. Four-liter reagent and waste bottles can crack and leak. Each 4-liter bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position. Wear appropriate eyewear, clothing, and gloves when handling reagent and waste bottles.



WARNING! CHEMICAL STORAGE HAZARD. Never collect or store waste in a glass container because of the risk of breaking or shattering. Reagent and waste bottles can crack and leak. Each waste bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position. Wear appropriate eyewear, clothing, and gloves when handling reagent and waste bottles.

Chemical safety guidelines

To minimize the hazards of chemicals:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. (See [“About SDSs” on page 138.](#))
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the SDS.
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the SDS.
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer’s cleanup procedures as recommended in the SDS.
- Comply with all local, state/provincial, or national laws and regulations related to chemical storage, handling, and disposal.



SDSs

About SDSs Chemical manufacturers supply current Safety Data Sheets (SDSs) with shipments of hazardous chemicals to new customers. They also provide SDSs with the first shipment of a hazardous chemical to a customer after an SDS has been updated. SDSs provide the safety information you need to store, handle, transport, and dispose of the chemicals safely.

Each time you receive a new SDS packaged with a hazardous chemical, be sure to replace the appropriate SDS in your files.

Obtaining SDSs The SDS for any chemical supplied by Applied Biosystems is available to you free 24 hours a day. To obtain SDSs:

1. Go to www.appliedbiosystems.com, click **Support**, then select **SDS**.
2. In the Keyword Search field, enter the chemical name, product name, SDS part number, or other information that appears in the SDS of interest. Select the language of your choice, then click **Search**.
3. Find the document of interest, right-click the document title, then select any of the following:
 - **Open** – To view the document
 - **Print Target** – To print the document
 - **Save Target As** – To download a PDF version of the document to a destination that you choose



Note: For the SDSs of chemicals not distributed by Applied Biosystems, contact the chemical manufacturer.



Chemical waste safety

Chemical waste hazards



CAUTION! HAZARDOUS WASTE. Refer to Safety Data Sheets and local regulations for handling and disposal.



WARNING! CHEMICAL WASTE HAZARD. Wastes produced by Applied Biosystems instruments are potentially hazardous and can cause injury, illness, or death.



WARNING! CHEMICAL STORAGE HAZARD. Never collect or store waste in a glass container because of the risk of breaking or shattering. Reagent and waste bottles can crack and leak. Each waste bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position. Wear appropriate eyewear, clothing, and gloves when handling reagent and waste bottles.

Chemical waste safety guidelines

To minimize the hazards of chemical waste:

- Read and understand the Safety Data Sheets (SDSs) provided by the manufacturers of the chemicals in the waste container before you store, handle, or dispose of chemical waste.
- Provide primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the SDS.
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the SDS.
- Handle chemical wastes in a fume hood.
- After emptying a waste container, seal it with the cap provided.
- Dispose of the contents of the waste tray and waste bottle in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations.

Waste disposal

If potentially hazardous waste is generated when you operate the instrument, you must:

- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure the health and safety of all personnel in your laboratory.



- Ensure that the instrument waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- ⓘ **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.

Biological hazard safety

General biohazard



WARNING! BIOHAZARD. Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Follow all applicable local, state/provincial, and/or national regulations. Wear appropriate protective equipment, which includes but is not limited to: protective eyewear, face shield, clothing/lab coat, and gloves. All work should be conducted in properly equipped facilities using the appropriate safety equipment (for example, physical containment devices). Individuals should be trained according to applicable regulatory and company/institution requirements before working with potentially infectious materials. Read and follow the applicable guidelines and/or regulatory requirements in the following:

- U.S. Department of Health and Human Services guidelines published in *Biosafety in Microbiological and Biomedical Laboratories* (stock no. 017-040-00547-4; www.cdc.gov/od/ohs/biosfty/bmbl5/bmbl5toc.htm)
- Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR§1910.1030; www.access.gpo.gov/nara/cfr/waisidx_01/29cfr1910a_01.html).
- Your company's/institution's Biosafety Program protocols for working with/handling potentially infectious materials.

Additional information about biohazard guidelines is available at:

www.cdc.gov





3'-end modification	Process by which dUTP is added to the 3' end of the P2 Adaptor on the templated beads using a terminal transferase reaction
aqueous phase	Emulsion component comprised of SOLiD™ P1 DNA Beads, library template, primers, DNA polymerase, dNTPs, and water
broken emulsion	An emulsion in which the aqueous phase appears at the bottom of a PCR plate well prior to emulsion break
Denaturing Buffer solution (or prepared Denaturing Buffer solution)	Solution made up of Denaturing Buffer and Denaturant used to make the template on templated beads single-stranded and to dissociate templated beads from enrichment beads during the enrichment step
emulsion break	Process by which the micro-reactors in an emulsion are broken using 2-butanol to allow processing of templated beads
emulsion PCR (ePCR)	Process by which DNA fragments are clonally amplified onto beads in individual droplets in an emulsion
enrichment	Process by which templated beads are isolated from non-amplifying beads using enrichment beads
enrichment beads	Polystyrene beads with a single-stranded P2 Adaptor attached to capture templated beads
full-scale templated bead preparation	Templated bead preparation process that yields 150 to 300 million templated beads
library	Set of DNA tags prepared from the same biological sample to be sequenced on the SOLiD™ System
macro-scale templated bead preparation	Templated bead preparation process that yields 600 million to 2.4 billion templated beads
micro-reactor	Droplet of aqueous phase in the emulsion in which amplification takes place
mini-scale templated bead preparation	Templated bead preparation process that yields 75 to 150 million templated beads
monoclonal bead	Templated bead with a single template

monoclonal micro-reactor	Micro-reactor containing a single template
monoclonal bead	Templated bead with a single template
multi-bead micro-reactor	Micro-reactor containing multiple beads
non-amplifying bead	SOLiD™ P1 DNA bead with no template
non-clonal micro-reactor	Micro-reactor containing no template
oil phase	Emulsion component of oil and emulsifiers
optimal library concentration	Library template concentration that gives the best sequencing results
P2-enriched beads	Enriched, templated beads
polyclonal bead	Templated bead with multiple templates
pulse-spin	Place the tube in a microfuge and spin for a few seconds to bring down any beads or liquid stuck on the walls of the tube.
remove the supernatant	Use a pipette to carefully remove the liquid from the tube without disturbing any beads.
resuspend the beads	<p>The beads can be resuspended in one of two ways:</p> <ul style="list-style-type: none"> • Gently pipet the solution up and down until the beads are suspended. Using a slower speed to aspirate and expel the solution minimizes the amount of beads that stick to the inside of the pipette tip. • Vortex the solution until all of the beads are suspended. Place the beads in a microfuge, then pulse-spin the beads for a few seconds to bring down any beads stuck on the walls of the tube. Do not over-spin the beads, or the beads pellet.
SOLiD™ P1 DNA Beads	Bead with P1 Adaptor attached
sonicate the beads	Place the tube containing the beads in the appropriate tube holder, then place the holder in the Covaris™ S2 System. Next, run the appropriate Covaris™ S2 program.
templated bead preparation	Process of adding library template to beads by emulsion PCR, enriching the beads to remove beads without template, and modifying the 3' end of the template on the beads to prepare for bead deposition and sequencing
templated beads	SOLiD™ P1 DNA Beads with amplified library template attached



titration	Library template concentration used to prepare an emulsion
workflow analysis (WFA) run	Type of run on the SOLiD™ System in which a small portion of templated beads are deposited and analyzed to test for templated bead quality



Related documentation

Document	Part number	Description
<i>Applied Biosystems SOLiD™ 4 System Library Preparation Guide</i>	4445673	Describes how to prepare libraries.
<i>Applied Biosystems SOLiD™ 4 System Library Preparation Quick Reference Card</i>	4445674	Provides brief, step-by-step procedures for preparing libraries.
<i>Applied Biosystems SOLiD™ 4 System Templated Bead Preparation Quick Reference Card</i>	4448329	Provides brief, step-by-step procedures for preparing templated beads by emulsion PCR (ePCR), required before sequencing on the SOLiD™ 4 System.
<i>Applied Biosystems SOLiD™ 4 System Instrument Operation Guide</i>	4448379	Describes how to load and run the SOLiD™ 4 System for sequencing.
<i>Applied Biosystems SOLiD™ 4 System Instrument Operation Quick Reference Card</i>	4448380	Provides brief, step-by-step procedures for loading and running the SOLiD™ 4 System.
<i>Applied Biosystems SOLiD™ 4 System Site Preparation Guide</i>	4448639	Provides all the information that you need to set up the SOLiD™ 4 System.
<i>Applied Biosystems SOLiD™ 4 System SETS Software User Guide</i>	4448411	Provides an alternate platform to monitor runs, modify settings and reanalyze previous runs that are performed on the SOLiD System.
<i>Applied Biosystems SOLiD™ 4 System ICS Software Help</i>	—	Describes the software and provides procedures for common tasks (see the Instrument Control Software).
<i>BioScope™ Software for Scientists Guide</i>	4448431	Provides a bioinformatics analysis framework for flexible application analysis (data-generated mapping, SNPs, count reads) from sequencing runs.
<i>Working with SOLiDBioScope.com™ Quick Reference Card</i>	4452359	Provides an online suite of software tools for Next Generation Sequencing (NGS) analysis. SOLiDBioScope.com™ leverages the scalable resources of cloud computing to perform compute-intensive NGS data processing.
<i>Applied Biosystems SOLiD™ 4 System Software Integrated Workflow Quick Reference Guide</i>	4448432	Describes the relationship between the softwares comprising the SOLiD 4 platform and provides quick step procedures on operating each software to perform data analysis.
<i>Applied Biosystems SOLiD™ 4 System Product Selection Guide</i>	4452360	Provides a quick guide to the sequencing kits you need to perform fragment, paired end, mate-pair, multiplex fragment, and multiplex paired end sequencing.

Document	Part number	Description
<i>Applied Biosystems SOLiD™ System SOLiD™ 3 Plus to SOLiD™ 4 System User Documentation Changes</i>	4451929	Provides a brief summary of changes made between the SOLiD™ 3 Plus System documentation and the SOLiD™ 4 System documentation.
<i>Applied Biosystems SOLiD™ 4 Upgrade Checklist</i>	4449773	Provides a checklist to ensure that all necessary preparations are made before upgrading to the SOLiD™ 4 System and provides a list of orderable consumables.



Note: For additional documentation, see [“How to obtain support”](#) on page 8.



Send us your comments

Applied Biosystems welcomes your comments and suggestions for improving its user documents. You can e-mail your comments to:

techpubs@appliedbiosystems.com

IMPORTANT! The e-mail address above is for submitting comments and suggestions relating *only* to documentation. To order documents, download PDF files, or for help with a technical question, see [“How to obtain support” on page 8](#).



Documentation
Send us your comments

A

Applied Biosystems
 customer feedback on documentation 149
 Information Development department 149

B

biohazardous waste, handling 141
 bold text, when to use 8

C

CAUTION, description 7
 checklists and workflow tracking forms 119
 chemical safety 137
 chemical waste safety 139
 conventions
 bold text 8
 for describing menu commands 8
 IMPORTANTs! 8
 in this guide 8
 italic text 8
 Notes 8
 user attention words 8
 Covaris™ S2 System 125
 customer feedback, on Applied Biosystems documents 149

D

DANGER, description 7
 documentation 147
 documentation, related 147

F

formulas and calculations 117

G

glossary 143
 guidelines
 chemical safety 137
 chemical waste disposal 139
 chemical waste safety 139

H

hazards. *See* safety

I

IMPORTANT, description 7
 Information Development department, contacting 149
 instrument operation, safety 136
 instrument warranty 129
 introduction 9
 italic text, when to use 8

M

macro-scale
 4 ePCR reactions 76
 8 ePCR reactions 86
 menu commands, conventions for describing 8
 MSDSs. *See* SDS

P

physical hazard safety 136
 preface 7
 prepare templated beads 21, 45, 69, 75, 85
 full-scale 46
 macro-scale 70
 mini-scale 22
 pressurized fluids, safety 136
 program the Eppendorf Repeater® Xstream Pipettor 106

Q

quantitate the beads using a hemocytometer 115
 quantitate the beads using the NanoDrop® ND-1000 Spectrophotometer 107, 112

R

radioactive waste, handling 140
 required materials 95

S

safety 135
 before operating the instrument 136
 biological hazards 141
 chemical 137
 chemical waste 139
 guidelines 137, 139
 instrument operation 136
 physical hazard 136

pressurized fluids. *See also* compressed gases, safety
solvents [136](#)

SDSs

about [7](#)
description [138](#)
obtaining [8](#), [138](#)

solvents, safety [136](#)

supplemental procedures [105](#)

generate a standard curve for calculating bead
concentration [112](#)

program the Eppendorf Repeater® Xstream
Pipettor [106](#)

quantitate the beads using a hemocytometer [115](#)

quantitate the beads using the NanoDrop® ND-1000
Spectrophotometer [107](#)

T

templated bead, overview [15](#)

text conventions [8](#)

training, information on [8](#)

U

user attention words, described [8](#)

W

WARNING, description [7](#)

warranty [129](#)

waste disposal, guidelines [139](#)

waste profiles, description [139](#)

Part Number 4448378 Rev. B 03/2010



Applied Biosystems

850 Lincoln Centre Drive | Foster City, CA 94404 USA
Phone 650.638.5800 | Toll Free 800.345.5224
www.appliedbiosystems.com

Technical Resources and Support

For the latest technical resources and support information
for all locations, please refer to our Web site at
www.appliedbiosystems.com/support