

# Ion ReproSeq™ PGS Kits

## USER GUIDE

### Single cell library preparation and sequencing on the Ion PGM™ System

**Catalog Numbers** A28888, A28889, A28890, A29142, A29143, A29144

**Publication Number** MAN0013762

**Revision** F.0

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Revision history:** Pub. No. MAN0013762

Revision	Date	Description
F.0	7 November 2016	<ul style="list-style-type: none"><li>Recommended input volume of 10 pM library pool used in the IA reaction increased from 10 <math>\mu</math>L to 18 <math>\mu</math>L.</li><li>Topic organization in Chapter 4, "Prepare template-positive Ion PGM™ Template IA ISPs" restructured for ease of use.</li></ul>
E.0	6 April 2016	<ul style="list-style-type: none"><li>Reagent cap colors in Ion PGM™ Template IA Reagents 500 and Reactions 500 updated.</li><li>Graphics enhanced.</li></ul>
D.0	12 January 2016	<ul style="list-style-type: none"><li>Maximum recommended libraries loaded on one Ion 316™ chip changed from 16 to 15.</li><li>Minor clarifications made in the Template IA workflow.</li><li>Clarifications made in the sequencing workflow, including use of the Line Clear feature in initialization.</li><li>Procedure added for editing the Ion Reporter™ Ion ReproSeq™ workflow with a different CNV Confidence Range.</li><li>Appendix on Ion OneTouch™ ES Instrument setup and maintenance deleted. Users referred to the <i>Ion OneTouch™ 2 System User Guide</i> (Pub. No. MAN0014388).</li></ul>
C.0	7 October 2015	<ul style="list-style-type: none"><li>Minor clarifications made in the Ion SingleSeq™ library preparation workflow.</li><li>Updated for Ion Reporter™ software v5.0, including:<ul style="list-style-type: none"><li>Ion Reporter™ analysis simplified with availability of pre-loaded Ion ReproSeq™ control baseline and aneuploidy workflow.</li><li>Instructions added for use of IGV Light in CNV analysis.</li></ul></li></ul>
B.0	9 August 2015	Configurations added for Ion ReproSeq™ PGS 314 and 316 kits, with and without chips.
A.0	17 July 2015	New User Guide. Provides detailed, step-by-step instructions on how to use the Ion ReproSeq™ PGS Kit to extract gDNA from single cells, prepare libraries, and generate templated Ion Sphere™ Particles for sequencing using the Ion PGM™ System.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

**Corporate entity:** Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

**Notice to purchaser:** The Ion SingleSeq™ Kit contains PicoPLEX™ technology developed and manufactured by Rubicon Genomics, Inc., Ann Arbor, Michigan, USA. PicoPLEX™ technology is covered by applicable US patents and corresponding patents outside the US.

**Trademarks:** All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Rainin and Pipet-Lite are trademarks of Rainin Instrument, LLC. ELGA and PURELAB are trademarks of VWS (UK) Ltd. iPod is a trademark of Apple, Inc. Agilent, Bioanalyzer, Mx3000P, Mx3005P, and Mx4000 are trademarks of Agilent Technologies, Inc. Agencourt and AMPure are trademarks of Beckman Coulter, Inc. Guava and easyCyte are trademarks of Guava Technologies, Inc. Kimwipes is a trademark of Kimberly-Clark Corporation. Eppendorf and Eppendorf LoBind are trademarks of Eppendorf AG. Luer-Lok is a trademark of Becton, Dickinson and Company. Milli-Q is a trademark of Millipore Corporation. Tween is a trademark of Uniquema Americas LLC. Xiameter is a trademark of Dow Corning Corporation. PicoPLEX is a trademark of Rubicon Genomics, Inc.

©2016 Thermo Fisher Scientific Inc. All rights reserved.

# Contents

About this guide .....	8
Purpose .....	8
Prerequisites .....	8
■ <b>CHAPTER 1</b> Product information .....	9
Product description .....	9
Software compatibility .....	9
Kit contents and storage .....	10
Kit configurations .....	10
Ion SingleSeq™ Kit .....	11
Ion PGM™ Template IA 500 Kit .....	13
Ion PGM™ Hi-Q™ Sequencing Kit .....	14
Ion PGM™ Wash 2 Bottle Kit .....	15
Ion Chip Kits .....	15
Required materials not supplied .....	16
Optional materials and equipment .....	18
Workflow overview .....	19
■ <b>CHAPTER 2</b> Before you begin .....	20
Before you begin .....	20
Update the software .....	20
Instrument installation by trained personnel only .....	20
Nucleic acid contamination .....	20
CO <sub>2</sub> contamination .....	20
Instrument vibration and clearances .....	21
Static electricity .....	21
Ventilation requirements .....	21
Gas cylinders .....	21
Perform a leak test .....	22
General procedural guidelines .....	23
Guidelines for the number of libraries per run .....	23
Guidelines for chip handling and use .....	23

■	<b>CHAPTER 3</b>	<b>Prepare and pool Ion SingleSeq™ libraries</b>	<b>24</b>
	Extract and amplify the gDNA		24
	Materials required		24
	Guidelines for sample handling		25
	Extract genomic DNA		25
	Pre-amplify the gDNA		26
	Amplify the libraries		27
	Library pooling, purification, and quantification		30
	Materials required		30
	Pool the libraries		30
	Purify the library pool		31
	Quantify the library pool		32
■	<b>CHAPTER 4</b>	<b>Prepare template-positive Ion PGM™ Template IA ISPs</b>	<b>34</b>
	Materials required		34
	Isothermal amplification (IA) reaction procedural guidelines		35
	Guidelines for preventing contamination		35
	IA reaction guidelines		35
	Before you begin		35
	Prepare the final library pool dilution		36
	Perform the IA reaction		36
	Recover the template-positive ISPs		38
■	<b>CHAPTER 5</b>	<b>Enrich the template-positive Ion PGM™ Template IA ISPs</b>	<b>39</b>
	Materials required		39
	Determine if a residual volume test is necessary		40
	Prepare reagents then fill the 8-well strip		41
	Prepare Melt-Off Solution		41
	Wash and resuspend the Dynabeads™ MyOne™ Streptavidin C1 Beads		41
	Fill the 8-well strip		42
	Prepare the Ion OneTouch™ ES		44
	Perform the Ion OneTouch™ ES run		46
	Sequence or store the template-positive ISPs		48
	Perform quality control on enriched ISPs		48

■	<b>CHAPTER 6</b>	<b>Create a Planned Run</b>	<b>49</b>
	About Planned Runs		49
	Create a Planned Run		50
	Planned Run wizard: key fields		53
■	<b>CHAPTER 7</b>	<b>Clean and initialize the Ion PGM™ Sequencer</b>	<b>54</b>
	Condition the Wash 2 Bottle for first use		54
	Cleaning		54
	Materials required		54
	Cleaning schedule		55
	Cleaning setup		55
	18 MΩ water cleaning		55
	Chlorite cleaning		56
	Initialize the Ion PGM™ System		57
	Materials required		57
	Initialization guidelines		58
	Before initialization		58
	Prepare the Wash 2 Bottle		58
	Prepare the Wash 1 and Wash 3 Bottles		60
	Begin the initialization		60
	Prepare the 50-mL Reagent Bottles with dNTPs		62
	Attach the sipper tubes and Reagent Bottles		63
■	<b>CHAPTER 8</b>	<b>Load the chip and start the sequencing run</b>	<b>64</b>
	Materials required		64
	Guidelines for sequencing		65
	Before you begin		65
	Optional: Prepare Ion Sphere™ Test Fragments		65
	Add controls to the enriched, template-positive ISPs		65
	Anneal the Sequencing Primer		66
	Perform Chip Check		67
	Bind the Sequencing Polymerase to the ISPs		69
	Prepare and load the chip		70
	Remove liquid from the chip		70
	Load the chip		71
	Select the Planned Run and perform the run		74
	Select the Planned Run		74
	Perform the run		75

■	<b>CHAPTER 9 Analyze the run</b>	<b>76</b>
	Launch an Ion Reporter™ analysis	76
	Edit the Ion Reporter™ ReproSeq Low-pass whole-genome aneuploidy workflow	78
	Visualize results with IGV Light viewer	81
■	<b>APPENDIX A Troubleshooting</b>	<b>84</b>
	Ion SingleSeq™ library preparation	84
	Ion OneTouch™ ES	84
	Chip Check	89
	Chip calibration (before loading sample)	90
	Chip calibration (after loading sample)	92
	Initialization—General errors	93
	Initialization—Auto pH errors	94
	Initialization—Reagent pH verification	99
	Ion Reporter™ results	100
■	<b>APPENDIX B Additional instrument information</b>	<b>101</b>
	Ion PGM™ Sequencer input and output connections	101
	Power the Ion PGM™ Sequencer on or off	102
	Power on	102
	Power off	102
	Update the Ion PGM™ System software	102
■	<b>APPENDIX C Manually adjust W2 pH</b>	<b>104</b>
	Materials and equipment needed	104
	Procedure	104
■	<b>APPENDIX D Safety</b>	<b>106</b>
	Symbols on this instrument	106
	Safety alerts on this instrument	107
	Safety information for instruments not manufactured by Thermo Fisher Scientific	108
	Instrument safety	108
	General	108
	Physical injury	108
	Electrical	108
	Cleaning and decontamination	109
	Laser	109

Safety and electromagnetic compatibility (EMC) standards .....	109
Safety .....	109
EMC .....	110
Environmental design .....	110
Chemical safety .....	111
Biological hazard safety .....	112
<b>Documentation and support .....</b>	<b>113</b>
Customer and technical support .....	113
Limited product warranty .....	113

# About this guide

---

**IMPORTANT!** Before using this product, read and understand the information in the “Safety” appendix in this document.

---

## Purpose

The *Ion ReproSeq™ PGS Kits User Guide* (Pub. No. MAN0013762) provides protocols and reference information for using an Ion ReproSeq™ PGS Kit to prepare whole genome pooled libraries from single cell samples, and prepare template-positive Ion Sphere™ Particles from them using an isothermal amplification reaction. A protocol for using the Ion PGM™ Hi-Q™ Sequencing Kit with Ion 314™, Ion 316™ and Ion 318™ Chips on the Ion PGM™ System is also provided in this guide.

## Prerequisites

This guide also assumes that you have:

- A general understanding of Ion Torrent™ sequencing chemistry and workflow
- Knowledge of techniques for handling and preparing DNA libraries





# Product information

## Product description

The Ion ReproSeq™ PGS Kits are bundles of four or five sub-kits which, when used in conjunction with the Ion PGM™ System, provide reagents and materials for whole genome amplification and sequencing to detect chromosomal aneuploidies, chromosome arm events (>48 Mb), and copy number variations starting from a single cell.

The Ion SingleSeq™ Kit (24 reactions/kit) include reagents to extract, amplify, and barcode genomic DNA. The Ion PGM™ Template IA 500 Kit includes reagents for preparing 4 reactions of template-positive Ion PGM™ Template IA Ion Sphere™ Particles (ISPs) for sequencing with the Ion PGM™ System. Template preparation is performed using Ion IA technology, whereby DNA is clonally amplified onto a bead surface through a non-emulsion, isothermal reaction. The Ion PGM™ Hi-Q™ Sequencing Kit, and Ion PGM™ Wash 2 Bottle Kit include reagents and materials for 4 initializations and up to 12 sequencing runs on the Ion PGM™ Sequencer.

## Software compatibility

Ion ReproSeq™ PGS Kits are compatible with Torrent Suite™ Software v5.0 and later. Be sure to update your Torrent Server to the latest available version of Torrent Suite™ Software before using these kits. For data analysis, Ion Reporter™ Software v5.2 or later is recommended.

## Kit contents and storage

### Kit configurations

Six configurations of the Ion ReproSeq™ PGS Kit are available. The kits vary in the number of Ion SingleSeq™ Kits provided, the maximum number of samples processed per kit, the type of Ion sequencing chip provided, and whether Ion sequencing chips are included.

<b>Ion ReproSeq™ PGS Kit</b>	<b>Cat. No.</b>	<b>Samples per kit</b>
Ion ReproSeq™ PGS 314 Kit, with chips <ul style="list-style-type: none"> <li>• Ion SingleSeq™ Kit (1 Kit)</li> <li>• Ion PGM™ Template IA 500 Kit</li> <li>• Ion PGM™ Hi-Q™ Sequencing Kit</li> <li>• Ion PGM™ Wash 2 Bottle Kit</li> <li>• Ion 314™ Chip Kit v2 BC (8 chips)</li> </ul>	A28888	16 <sup>[1]</sup>
Ion ReproSeq™ PGS 316 Kit, with chips <ul style="list-style-type: none"> <li>• Ion SingleSeq™ Kit (5 kits)</li> <li>• Ion PGM™ Template IA 500 Kit</li> <li>• Ion PGM™ Hi-Q™ Sequencing Kit</li> <li>• Ion PGM™ Wash 2 Bottle Kit</li> <li>• Ion 316™ Chip Kit v2 BC (8 chips)</li> </ul>	A28889	120
Ion ReproSeq™ PGS 318 Kit, with chips <ul style="list-style-type: none"> <li>• Ion SingleSeq™ Kit (8 kits)</li> <li>• Ion PGM™ Template IA 500 Kit</li> <li>• Ion PGM™ Hi-Q™ Sequencing Kit</li> <li>• Ion PGM™ Wash 2 Bottle Kit</li> <li>• Ion 318™ Chip Kit v2 BC (8 chips)</li> </ul>	A28890	192
Ion ReproSeq™ PGS 314 Kit, without chips <ul style="list-style-type: none"> <li>• Ion SingleSeq™ Kit (1 Kit)</li> <li>• Ion PGM™ Template IA 500 Kit</li> <li>• Ion PGM™ Hi-Q™ Sequencing Kit</li> <li>• Ion PGM™ Wash 2 Bottle Kit</li> </ul>	A29142	16
Ion ReproSeq™ PGS 316 Kit, without chips <ul style="list-style-type: none"> <li>• Ion SingleSeq™ Kit (5 kits)</li> <li>• Ion PGM™ Template IA 500 Kit</li> <li>• Ion PGM™ Hi-Q™ Sequencing Kit</li> <li>• Ion PGM™ Wash 2 Bottle Kit</li> </ul>	A29143	120

Ion ReproSeq™ PGS Kit	Cat. No.	Samples per kit
Ion ReproSeq™ PGS 318 Kit, without chips <ul style="list-style-type: none"> <li>• Ion SingleSeq™ Kit (8 kits)</li> <li>• Ion PGM™ Template IA 500 Kit</li> <li>• Ion PGM™ Hi-Q™ Sequencing Kit</li> <li>• Ion PGM™ Wash 2 Bottle Kit</li> </ul>	A29144	192

<sup>[1]</sup> Although an Ion SingleSeq™ kit provides reagents for preparing 24 barcoded libraries, we recommend that no more than 2 libraries be loaded on a Ion 314™ Chip v2. See "Guidelines for the number of libraries per run" on page 23 for the recommended maximum number of libraries per sequencing run for each kind of chip.

## Ion SingleSeq™ Kit

Use the Ion SingleSeq™ Kit (Part No. A28955; 24 reactions/kit) to extract, amplify, and barcode genomic DNA. Immediately before use, thaw non-enzyme tubes on ice as needed.

Component	Amount	Storage
Cell Extraction Buffer (green cap)	120 µL	-30°C to -10°C
Extraction Enzyme Dilution Buffer (violet cap)	115.2 µL	
Cell Extraction Enzyme (yellow cap)	4.8 µL	
Pre-Amplification Buffer (red cap)	115.2 µL	
Pre-Amplification Enzyme (white cap)	4.8 µL	
Amplification Buffer (orang cap)	648 µL	
Amplification Enzyme (blue cap)	12 µL	
Nuclease-free Water (clear cap)	108 µL	
Human CEPH Genomic DNA Control, 100 µg/mL <sup>[1]</sup> (red cap)	30 µL	
Barcode Plate (Ion SingleSeq™ Barcode Set 1 loaded in wells A1–H3)	1 plate; 5 µL/well	

<sup>[1]</sup> Requires dilution prior to use in troubleshooting.

**About the Ion SingleSeq™ Barcode Set 1**

The Ion SingleSeq™ Kit is designed for high throughput applications and includes a 96-well Barcode Plate containing 24 unique Barcode Adapters. Each well contains sufficient volume for a single use. The plate is sealed with foil that can be pierced with a pipet tip to collect the required amount of the Barcode Adapter for a reaction.

To minimize cross-contamination, either use all 24 barcodes at once, or if you are using a subset, seal the pierced wells with laboratory tape. Do not thaw and re-freeze the plate more than 4 times.

**IMPORTANT!**

- Do not substitute barcode oligonucleotides from any other source for the Barcode Adapters contained in this kit.
- Avoid repeated freezing and thawing of the plate. Do not exceed 4 freeze/thaw cycles.

	1	2	3	4	5	6	7	8	9	10	11	12
A	1	9	17									
B	2	10	18									
C	3	11	19									
D	4	12	20									
E	5	13	21									
F	6	14	22									
G	7	15	23									
H	8	16	24									

Barcode Plate containing the Ion SingleSeq™ Barcode Set 1. Twenty-four Barcode Adapters are loaded in wells A1 through H3.

## Ion PGM™ Template IA 500 Kit

**IMPORTANT!** Use only the Ion PGM™ Template IA 500 Kit with the procedure in this guide. Do not use the Ion OneTouch™ 2 Instrument. Do not mix reactions or disposables, including plates, solutions, and kit reagents from other template preparation kits. We have verified this protocol using this specific material. Substitution can adversely affect performance.

**IMPORTANT!** Ion PGM™ Template IA Reactions 500 is shipped at 4°C to 8°C. On receipt, store at –30°C to –10°C. Immediately before use, thaw tubes on ice as needed.

### Kit contents

Contents <sup>[1]</sup>	Amount	Storage
Ion PGM™ Template IA Supplies 500 (Part No. A24618)		
Ion OneTouch™ ES Supplies, including: <ul style="list-style-type: none"><li>Eppendorf™ LoRetention Dualfilter, 300 µL PCR pipette tips (5)</li><li>8-well strips (12)</li></ul>	1 bag	15°C to 30°C
Ion PGM™ Template IA Reagents 500 (Part No. A24619)		
Ion PGM™ Template IA Pellets 500	4 foil pouches <sup>[2]</sup>	4°C to 8°C
Ion PGM™ Template IA ISP Dilution Buffer (yellow cap)	1 mL	
Ion PGM™ Template IA Start Solution (purple cap)	2 × 1200 µL	
Ion PGM™ Template IA Reactions 500 (Part No. A24620)		
Ion PGM™ Template IA Ion Sphere™ Particles (orange cap)	86 µL	–30°C to –10°C
Ion PGM™ Template IA Primer Mix S (black cap)	40 µL	
Ion PGM™ Template IA Primer Mix L (blue cap)	40 µL	
Ion PGM™ Template IA Rehydration Buffer (white cap)	2 × 1.5 mL	
Ion PGM™ Template IA Solutions 500 (Part No. A24621)		
Ion PGM™ Template IA Stop Solution	2.8 mL	15°C to 30°C
Ion PGM™ Template IA Recovery Solution	7 mL	
Ion PGM™ Template IA Wash Solution	4 mL	
MyOne™ Beads Wash Solution (green cap)	1.04 mL	
Neutralization Solution (red cap)	40 µL	
Tween™ Solution	3.5 mL	

<sup>[1]</sup> We have verified this protocol using this specific material. Substitution can adversely affect performance.

<sup>[2]</sup> One Ion PGM™ Template IA Pellet 500 per tube per pouch. One pellet used per IA 500 amplification reaction.

**Ion PGM™ Hi-Q™  
Sequencing Kit**

The Ion PGM™ Hi-Q™ Sequencing Kit (Cat. No. A25592) contains sufficient reagents and materials for 4 initializations of the Ion PGM™ Sequencer. Using 250 flows per run, the kit enables up to 3 runs per initialization, and up to 12 runs per kit.<sup>[1]</sup> The kit includes the following components:

Component	Amount	Storage
Ion PGM™ Sequencing Supplies (Part No. A25587)		
Wash 1 Bottle w/ label (250 mL)	1 bottle	15°C to 30°C
Wash 3 Bottle w/ label (250 mL)	1 bottle	
Ion PGM™ Reagent Bottle Sipper Tubes (blue)	16 tubes	
Ion PGM™ Wash Bottle Sipper Tubes (gray)	8 tubes for 250-mL bottles 4 tubes for 2-L bottles	
Reagent Bottles w/ labels (50 mL)	25 bottles	
Ion PGM™ Hi-Q™ Sequencing Reagents (Part No. A25588)		
Ion PGM™ Hi-Q™ Sequencing Polymerase (yellow cap)	36 µL	-30°C to -10°C
Sequencing Primer (white cap)	144 µL	
Control Ion Sphere™ Particles (clear cap)	60 µL	
Ion PGM™ Hi-Q™ Sequencing Solutions (Part No. A25589)		
Ion PGM™ Hi-Q™ Sequencing W2 Solution	4 × 126.25 mL	2°C to 8°C (store W2 Solution protected from light)
Ion PGM™ Cleaning Tablet	4 tablets	
Annealing Buffer	12 mL	
Ion PGM™ Hi-Q™ Sequencing W3 Solution	2 × 100 mL	
Ion PGM™ Hi-Q™ Sequencing dNTPs (Part No. A25590)		
Ion PGM™ Hi-Q™ Sequencing dGTP (black cap)	80 µL	-30°C to -10°C
Ion PGM™ Hi-Q™ Sequencing dCTP (blue cap)	80 µL	
Ion PGM™ Hi-Q™ Sequencing dATP (green cap)	80 µL	
Ion PGM™ Hi-Q™ Sequencing dTTP (red cap)	80 µL	

<sup>[1]</sup> If you are using a kit that provides chips, and you wish to perform 3 runs per initialization, you will need to order additional sequencing chips separately.

## Ion PGM™ Wash 2 Bottle Kit

The Ion PGM™ Wash 2 Bottle Kit (Cat. No. A25591) includes the following components:

Component	Amount	Storage
Wash 2 Bottle w/ label (2 L) <b>Note:</b> Must be conditioned at least 8 hours before use as described in "Condition the Wash 2 Bottle for first use" on page 54	1 bottle	15°C to 30°C
Wash 2 Bottle Conditioning Solution	125 mL	

## Ion Chip Kits

The following Ion Chip Kits are compatible with the Ion ReproSeq™ PGS Kits, and can be ordered separately if your Ion ReproSeq™ PGS Kit does not include chips:

Component	Quantity	Catalog No.	Storage
Ion 314™ Chip v2 BC	8 pack	4488144	15°C to 30°C
Ion 316™ Chip v2 BC	4 pack	4488145	
	8 pack	4488149	
Ion 318™ Chip v2 BC	4 pack	4488146	
	8 pack	4488150	

## Required materials not supplied

Unless otherwise indicated, all materials are available through **thermofisher.com**.  
MLS: Fisher Scientific (**fisherscientific.com**) or other major laboratory supplier.

Materials and equipment required for Ion SingleSeq™ library preparation.

✓	Item	Source
	ViiA™ 7 Real-Time PCR System with 96-Well Block (recommended), <i>or</i>	4453534
	Veriti™ 96-Well Thermal Cycler (0.2-mL block; for standard PCR amplification), <i>or</i>	4452300
	equivalent real-time or standard thermal cycler with heated lid	
	MicroAmp™ 96-Well Optical Reaction Plate	N8010560
	MicroAmp™ Optical Adhesive Film (for real-time PCR)	4360954
	MicroAmp™ Clear Adhesive Film (for standard endpoint PCR)	4306311
	<i>or equivalent</i>	
	<i>Optional:</i> MicroAmp™ Optical 8-Tube Strip; 0.2-mL <sup>[1]</sup>	4316567
	MicroAmp™ Optical 8-Cap Strip	4323032
	<i>or equivalent</i>	
	96-Well Tray/Retainer Set specific to your instrument	
	Agencourt™ AMPure™ XP Reagent	Beckman Coulter A63880
	DynaMag™ -2 magnet	12321D
	1.5-mL Eppendorf LoBind™ Tubes (DNA) <sup>[2]</sup>	Fisher Scientific 13-6987-91
	Qubit™ 3.0 Fluorometer <i>or</i>	Q33216
	Qubit™ 2.0 Fluorometer <sup>[3]</sup>	
	Qubit™ dsDNA HS Assay Kit	Q32851 Q32854
	Qubit™ Assay Tubes	Q32856
	Ethanol	MLS
	Low TE (10 mM Tris pH 8.0, 0.1 mM EDTA)	MLS
	Phosphate-buffered saline (1X PBS, Ca <sup>2+</sup> -free, Mg <sup>2+</sup> -free, BSA-free)	MLS
	Nuclease-free Water	AM9932
	SYBR™ Green I Nucleic Acid Gel Stain, 10,000X concentrate in DMSO <sup>[4]</sup>	S-7567
	ROX Reference Dye <sup>[4]</sup>	12223-012



- [1] For standard end-point PCR amplification, standard tube strip and caps may be substituted.  
 [2] May substitute 2.0-mL Eppendorf LoBind™ Tubes for 1.5-mL tubes.  
 [3] Supported but no longer available for purchase.  
 [4] Required if monitoring amplification by real-time PCR.

Materials and equipment required for templating.

✓	Item	Source
	Ion OneTouch™ ES Instrument	4473574
	2.0-mL Eppendorf LoBind™ Tubes (DNA)	Fisher Scientific 13-698-792
	Heat block set to 40°C	MLS
	Ion PGM™ Enrichment Beads (Dynabeads™ MyOne™ Streptavidin C1 Beads)	4478525 65001, 65002
	DynaMag™ -2 magnet	12321D

Materials and equipment required for sequencing.

✓	Item	Source
	Tank of compressed nitrogen (grade 4.5, 99.995% or better) (see "Gas cylinders" on page 21)	MLS
	Multistage (dual-stage) gas regulator (0-50 PSI, 2-3 Bar output)	Fisher Scientific NC0393866 or MLS
	<b>Choose from one of the following systems:</b> ELGA™ PURELAB™ Flex 3 Water Purification System Equivalent 18-MΩ water purification system	4474524 MLS
	Microcentrifuge (capable of >15,500 × g, fits 1.5-mL and 0.2-mL microcentrifuge tubes)	MLS
	0.22-μm or 0.45-μm vacuum filtration system and filters (nylon or PVDF filters, 1-L volume)	MLS
	Rainin™ Pipet-Lite™ LTS Pipette L-100XLS 10–100 μL <sup>[1]</sup> (Alternatives from Gilson and Eppendorf may be used)	Rainin 17014384
	Rainin™ Pipet-Lite™ LTS Pipette L-20XLS 2–20 μL <sup>[1]</sup> (Alternatives from Gilson and Eppendorf may be used)	Rainin 17014392
	Rainin™ LTS pipette tips, 200 μL, SR-L200F (Alternatives from Gilson and Eppendorf may be used)	Rainin 17005859
	Rainin™ LTS pipette tips, 20 μL, SR-L10F <sup>[1]</sup> (Alternatives from Gilson and Eppendorf may be used)	Rainin 17005860
	PCR tubes, Flat Cap, 0.2-mL (do not use polystyrene tubes)	Fisher Scientific 14-222-262

✓	Item	Source
	Vortexer with a rubber platform	MLS
	Thermal cycler with a heated lid	MLS
	Graduated cylinders (1 L or 2 L volume)	MLS
	Glass bottle (1 L)	MLS
	15-mL conical tubes	MLS
	NaOH (10 M), molecular biology grade	MLS
	Pipette set and filtered tips, P2, P20, P200, and P1000 µL	MLS
	Microcentrifuge tubes, 1.5-mL or 1.7-mL	MLS
	Syringe, 10 CC, Female Luer-Lok™ (used for clearing lines)	Provided with the Ion PGM™ Sequencer, or MLS
	Ion PGM™ 2.5 L Waste Bottle <sup>[2]</sup>	4482565
<b>Optional materials</b>		
	Ion PGM™ Controls Kit <sup>[3]</sup>	4480449
	Ion PGM™ Sequencing Sippers Kit <sup>[4]</sup>	4478682
	Non-interruptible Power Supply (UPS) <sup>[5]</sup>	MLS

<sup>[1]</sup> Ensure tips from any vendors are low binding tips. Required for loading samples onto the Ion PGM™ Chips.

<sup>[2]</sup> Required one-time purchase for older ("Orange") versions of the Ion PGM™ System.

<sup>[3]</sup> Not commonly needed, but available for troubleshooting.

<sup>[4]</sup> Contains additional sipper tubes; not commonly needed.

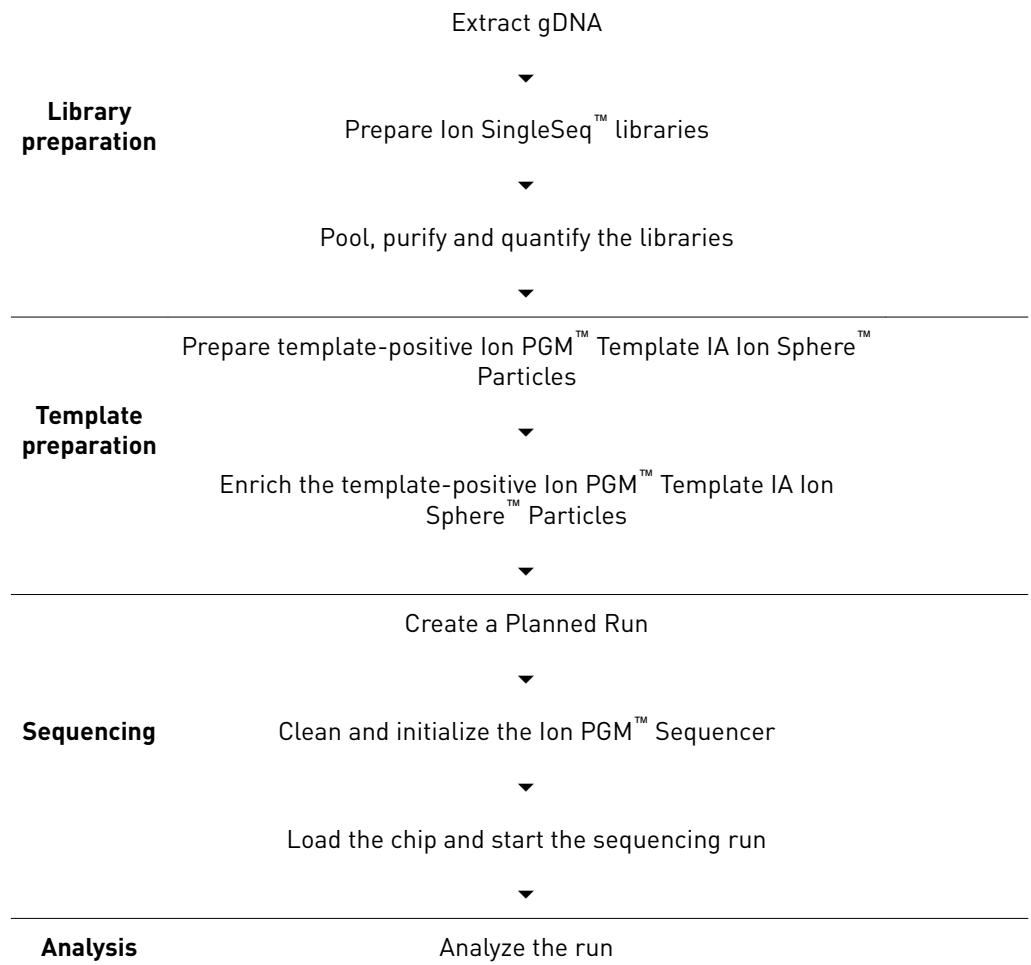
<sup>[5]</sup> We recommend using a non-interruptible power supply (UPS) for laboratories that experience frequent power outages or line voltage fluctuations. The UPS must be compatible with 1500 W output or higher. The 1500 VA unit from APC provides ~11 minutes of backup power for an Ion PGM™ System.

## Optional materials and equipment

The materials in the following table are optional and required only to verify and adjust the pH of the W2 Solution prior to sequencing the Ion chips.

✓	Item	Source
	Orion™ 3-Star Plus Benchtop pH Meter Kit with electrode, electrode stand, and calibration buffers (or equivalent)	Fisher Scientific 1112003
	1 N HCl	MLS
	Magnetic stirrer (must hold 2-L bottle)	MLS
	Magnetic stir bar (4 cm)	MLS
	Squirt bottle	MLS

## Workflow overview





# Before you begin

■ Before you begin .....	20
■ Gas cylinders .....	21
■ General procedural guidelines .....	23
■ Guidelines for chip handling and use .....	23

## Before you begin

For additional safety information, see Appendix D, “Safety”.

### Update the software

---

**IMPORTANT!** Before proceeding, make sure that you have updated the Torrent Suite™ and Ion PGM™ System software to Version 5.0 or later. See “Update the Ion PGM™ System software” on page 102.

---

### Instrument installation by trained personnel only

---

**IMPORTANT!** The Ion PGM™ System is installed by trained service personnel and must not be relocated without assistance from trained service personnel. See “Customer and technical support” on page 113.

---

### Nucleic acid contamination

---

**IMPORTANT!** A primary source of contamination is DNA fragments from previously processed samples. Do not introduce amplified DNA into the library preparation laboratory or work area.

---

---

**IMPORTANT!** Possible contamination can occur during the transfer of dNTPs into Reagent Tubes. Be careful to avoid cross contamination of dNTP stocks. Barrier tips are required for all pipetting steps. Change gloves after handling concentrated dNTP stocks.

---

### CO<sub>2</sub> contamination

---

**IMPORTANT!** Dry ice (solid CO<sub>2</sub>) must be kept away from areas where buffers, wash solutions, or sources of molecular biology grade water for the Ion PGM™ System are used. High air concentrations of subliming CO<sub>2</sub> may change the pH of such buffers during or after their preparation. The stability of the pH of these buffers is a critical factor in the performance of the Ion PGM™ System.

---

## Instrument vibration and clearances

---

**IMPORTANT!** Significant vibration during sequencing may add noise and reduce the quality of the measurements. The Ion PGM™ System must be installed on a bench that is free from vibrations or in contact with equipment that can cause vibrations to the bench (freezers, pumps, and other similar equipment).

---

---

**IMPORTANT!** Position the Ion PGM™ System so that the front bezel is a minimum of 12 in. (30.5 cm) and the Reagent Tubes containing dNTPs are a minimum of 8 in. (20.3 cm) from the front of the laboratory bench. Place the instrument at least 40 in. (1 meter) away from major sources of electronic noise such as refrigerators or microwaves.

---

## Static electricity

---

**IMPORTANT!** To avoid possible damage to the chip from static electricity, prior to handling chips, you must ground yourself on the grounding plate (located next to the chip clamp) by touching the grounding plate with your bare hand.

Do not place chips on non-grounded surfaces such as a bench.

Always use the grounding plate to hold chips that are not in the package inserted in the chip clamp or the Ion Chip™ Minifuge bucket.

---

## Ventilation requirements



**WARNING!** Instrumentation must be installed and operated in a well-ventilated environment, defined as having a minimum airflow of 6–10 air changes per hour. Assess the need for ventilation or atmospheric monitoring to avoid asphyxiation accidents from inert gases and/or oxygen depletion, and take measures to clearly identify potentially hazardous areas through training or signage. Please contact your Environmental Health and Safety Coordinator to confirm that the instruments will be installed and operated in an environment with sufficient ventilation.

---

## Gas cylinders

You must supply the required nitrogen gas cylinder and accessories for the installation. This instrument requires a pressurized house line or one size 1-A nitrogen gas cylinder that holds approximately 7.2 m<sup>3</sup> (257 ft<sup>3</sup>) of gas when full. Use only prepurified nitrogen of 99.995% (grade 4.5) or greater purity.



**CAUTION!** Damage to the instrument and its products can result from using impure gas, gases other than nitrogen, or an inadequate amount of gas.

---



**WARNING! EXPLOSION HAZARD.** Pressurized gas cylinders are potentially explosive. Always cap the gas cylinder when it is not in use, and attach it firmly to the wall or gas cylinder cart with approved brackets or chains.

---



**WARNING!** Gas cylinders are heavy and may topple over, potentially causing personal injury and tank damage. Cylinders should be firmly secured to a wall or work surface. Please contact your Environmental Health and Safety Coordinator for guidance on the proper installation of a gas cylinder.

---

## Perform a leak test

To perform a leak test on the gas cylinder:

1. Open the main tank shutoff valve. The high-pressure gauge of the gas tank regulator reads approximately 2,000–2,500 psi for a full tank.
2. Adjust the pressure to the instrument by slowly turning the pressure adjustment valve clockwise until the low-pressure gauge reads 30 psi.
3. Close the needle valve, then close the main tank valve.
4. Monitor the high-pressure gauge of the gas tank regulator for 5 minutes. There should be no noticeable drop in pressure.

If the pressure	Action
Drops in 5 minutes	There can be a leak at either the needle valve or the gas tank regulator itself. Check the fittings and resolve any problems, then continue with step 5.
Does not drop in 5 minutes	The instrument passes the leak test. Reopen the main tank valve and skip the following steps.

5. Open the main tank valve and the needle valve for at least 15 seconds to pressurize the instrument.
6. Close the main tank valve.
7. Monitor the high-pressure gas tank regulator gauge. There should be no more than a 100-psi drop in pressure after 5 minutes. Locate, then resolve any leaks. Turn the main tank valve back on.

## General procedural guidelines

- Use good laboratory practices to minimize cross-contamination of products. When designing the laboratory layout, consider the need for space separation of pre- and post-amplification activities. Dedicate laboratory supplies and/or equipment to the appropriate space to reduce significantly the potential for contamination.
- Pipet viscous solutions slowly and ensure complete mixing.
- Ensure that all reagents are completely thawed at room temperature, that is, no ice crystals are visible.
- Vortex all reagents, *except* for enzymes, for 5 seconds (mix enzymes by flicking the tube with your finger four times). Pulse centrifuge in a microcentrifuge for 3–5 seconds before use.

### Guidelines for the number of libraries per run

The maximum number of Ion SingleSeq™ libraries you can pool and sequence on one chip depends on the type of chip you use. Follow the guidelines in the table below:

Ion Chip	Recommended maximum libraries per run
Ion 314™ Chip v2 BC	2
Ion 316™ Chip v2 BC	15
Ion 318™ Chip v2 BC	24

## Guidelines for chip handling and use

- Always remove gloves before transferring chips onto or off the instrument. Hold chips by their edges when handling.
- To avoid damage due to electrostatic discharge (ESD), **do not place chips directly on the bench or any other surface**. Always place chips either on the grounding plate on the sequencer or in the Ion Chip™ Minifuge bucket.
- Used chips cannot be reused for sequencing. Used chips must be discarded or clearly marked for cleaning and initialization.



# Prepare and pool Ion SingleSeq<sup>TM</sup> libraries

■ Extract and amplify the gDNA .....	24
■ Library pooling, purification, and quantification .....	30

## Extract and amplify the gDNA

### Materials required    Provided in the Ion SingleSeq<sup>TM</sup> Kit (Part No. A28955):

- Cell Extraction Buffer
- Extraction Enzyme Dilution Buffer
- Cell Extraction Enzyme
- Pre-Amplification Buffer
- Pre-Amplification Enzyme
- Amplification Buffer
- Amplification Enzyme
- Nuclease-free Water
- Barcode Plate containing Ion SingleSeq<sup>TM</sup> Barcode Set 1

### Other materials:

- MicroAmp<sup>TM</sup> 96-well Optical Reaction Plate and adhesive film (or 0.2-mL 8-tube PCR tube strip with caps)
- Pipettor
- Pipette tips
- Real-time or standard PCR thermal cycler equipped with heated lid
- SYBR<sup>TM</sup> Green I and ROX dyes (Cat. Nos. S-7567 and 12223-012), if monitoring amplification by real-time PCR
- If performing troubleshooting, one of the following for diluting gDNA:
  - Low TE
  - PBS (Ca<sup>2+</sup>-, Mg<sup>2+</sup>-, BSA-free)
  - Nuclease-free Water



## Guidelines for sample handling

- **Keep the cells on ice.**
- For single cells or low-concentration input gDNA, do not insert pipette tip into sample. The single cell could adhere to the tip and get removed from the sample tube.
- Add liquids above the top of liquid in a tube, do not submerge the tip.
- **Do not vortex the cells.**
- After adding Extraction Enzyme Master Mix to samples, do not mix. Vortexing or pipetting up and down can cause loss of cells on the wall of the well or pipette tip.
- Keep cell lysis and amplification reactions on ice or a cold block during reaction setup, and keep Ion SingleSeq™ libraries on ice during library pooling and quantification.

## Extract genomic DNA

1. Prepare, by FACS or micro-manipulation, 1–10 cells per sample in up to 2.5 µL 1X PBS or Low TE, then pipet the cells into wells of a 96-well optical reaction plate.

**Note:** Instead of a 96-well reaction plate, a PCR tube strip can be used.

2. Add Cell Extraction Buffer (green cap) to each sample well to bring the total volume to 5 µL.

**Note:** You can also sort cells by FACS directly into Cell Extraction Buffer at a density of 1–10 cells/5 µL, then add 5 µL aliquots to reaction plate wells.

**Note:** If using gDNA for troubleshooting, mix 2 µL Human CEPH gDNA (red cap) at a concentration of 7.5 pg/µL with 3 µL Cell Extraction Buffer.

To dilute the 100 µg/mL gDNA stock solution to 7.5 pg/µL:

1. Add 2 µL of stock solution to 198 µL of Low TE. Vortex to mix, then centrifuge briefly.
2. Add 2 µL of this solution to 265 µL of Low TE. Vortex to mix, then centrifuge briefly.

Instead of Low TE, 1X PBS or Nuclease-free Water can be used.

3. (Optional) Prepare a Non-Template Control (NTC) by adding 2.5 µL 1X PBS to 2.5 µL Cell Extraction Buffer.
4. Prepare an Extraction Enzyme Master Mix in a 1.5-mL tube on ice according to the following table. Scale the volume according to the number of samples (N) and NTCs you have. Vortex briefly, then centrifuge to collect liquid at the bottom of the tube.

Extraction Enzyme Master Mix

Component	Volume per reaction	Volume for N reactions <sup>[1]</sup>
Extraction Enzyme Dilution Buffer (violet cap)	4.8 µL	$N \times 4.8 \mu\text{L} \times 1.1$
Cell Extraction Enzyme (yellow cap)	0.2 µL	$N \times 0.2 \mu\text{L} \times 1.1$

<sup>[1]</sup> 10% overage added to compensate for pipetting loss. Round to nearest 0.1 µL.

5. Add 5 µL Extraction Enzyme Master Mix to each cell or gDNA sample (10-µL final volume).

**Note:** Do not mix. Vortexing or pipetting up and down can cause loss of cells on the wall of the well or pipette tip.

6. Seal the plate, then centrifuge at  $1,000 \times g$  for 30 seconds to collect liquid at the bottom of the wells.
7. Incubate the samples in a thermal cycler using the following temperature program:

Temperature	Time
75°C	10 minutes
95°C	4 minutes
22°C	Hold

8. Remove the plate, centrifuge at  $1,000 \times g$  for 30 seconds to collect liquid at the bottom of the wells, and place on ice or a cold block.

Proceed to "Pre-amplify the gDNA".

## Pre-amplify the gDNA

1. Prepare a Pre-amplification Master Mix in a 1.5-mL tube on ice according to the following table. Scale the volume according to the number of samples (N) and NTCs you have. Vortex briefly and centrifuge for 30 seconds to collect liquid at the bottom of tube.

Pre-amplification Master Mix

Component	Volume per reaction	Volume for N reactions <sup>[1]</sup>
Pre-Amplification Buffer (red cap)	4.8 µL	$N \times 4.8 \mu\text{L} \times 1.1$
Pre-Amplification Enzyme (white cap)	0.2 µL	$N \times 0.2 \mu\text{L} \times 1.1$

<sup>[1]</sup> 10% overage added to compensate for pipetting loss. Round to nearest 0.1 µL.

2. Carefully remove the film seal from the plate, then add 5-µL Pre-amplification Master Mix to each sample well (15-µL final volume).  
**Note:** Do not mix. Vortexing or pipetting up and down can cause loss of DNA on the wall of the tube or pipette tip.
3. Apply a new adhesive film, then centrifuge at  $1,000 \times g$  for 30 seconds to collect liquid at the bottom of the wells.

4. Cycle samples in a thermal cycler according to the following program:

Step	Temperature	Time <sup>[1]</sup>	Number of cycles
1	95°C	2 minutes	1
2	95°C	15 seconds	12
	15°C	50 seconds	
	25°C	40 seconds	
	35°C	30 seconds	
	65°C	40 seconds	
	75°C	40 seconds	
3	4°C	Hold	1

<sup>[1]</sup> Cycling time is approximately 1 hour.

**IMPORTANT!** The thermal cycler must be equipped with a heated lid.

5. Remove the plate, centrifuge at  $1,000 \times g$  for 30 seconds to collect liquid at the bottom of the wells, and place on ice or a cold block.

Proceed to "Amplify the libraries".

## Amplify the libraries

You can perform the amplification reaction on a real-time PCR instrument to monitor amplification during the run. Monitoring can give an early indication of amplification failure in a particular sample, or that a library will be over- or under-represented after pooling.

1. If performing real-time PCR, prepare SYBR™ Green I/ROX dye mix:
  - a. Dilute the stock SYBR™ Green I reagent 1:1000 with Low TE buffer to make a 10X working solution.
  - b. Combine 10X SYBR™ Green I working solution and 25 µM ROX Reference Dye for the number of reactions (N) according to the following table:

Component	Volume per reaction	Volume for N reactions <sup>[1]</sup>
10X SYBR™ Green I working solution	0.5 µL	$N \times 0.5 \mu\text{L} \times 1.1$
25 µM ROX Reference Dye (or 2.5 µM, see below)	1.0 µL	$N \times 1.0 \mu\text{L} \times 1.1$
Nuclease-free Water (clear cap)	1.0 µL	$N \times 1.0 \mu\text{L} \times 1.1$

<sup>[1]</sup> 10% overage added to compensate for pipetting loss. Round to nearest 0.1 µL.

**IMPORTANT!** Use ROX Reference Dye at 25 µM for the following real-time PCR instruments: Applied Biosystems™ 7300, 7900HT, StepOne™, StepOnePlus™, ABI PRISM™ 7000, and 7700 instruments.

Dilute ROX Reference Dye to 2.5 µM with Nuclease-free Water and use 1.0 µL per reaction for the following instruments: Applied Biosystems™ 7500 and ViiA™ 7 instruments, and Agilent™ Mx3000P™, Mx3005P™, and Mx4000™ instruments.

**Note:** If performing standard endpoint PCR, use Nuclease-free Water instead of the dye mix, as indicated in the table in Step 3.

2. Prepare the Ion SingleSeq™ Barcode Set 1 plate:
  - a. Thaw the plate for 10 minutes at room temperature.
  - b. Centrifuge the plate briefly to collect contents at the bottom of the wells.
  - c. Wipe the foil seal with 70% ethanol, then allow it to dry.
3. Prepare an Amplification Master Mix in a 1.5-mL tube on ice according to the following table. Scale the volume according to the number (N) of samples and NTCs you have. Vortex briefly, then centrifuge to collect liquid at the bottom of tube.

Amplification Master Mix

Component	Volume per reaction	Volume for N reactions <sup>[1]</sup>
Amplification Buffer (orange cap)	27 µL	$N \times 27 \mu\text{L} \times 1.1$
Amplification Enzyme (blue cap)	0.5 µL	$N \times 0.5 \mu\text{L} \times 1.1$
SYBR™ Green I/ROX dye mix <sup>[2]</sup>	2.5 µL	$N \times 2.5 \mu\text{L} \times 1.1$

<sup>[1]</sup> 10% overage added to compensate for pipetting loss. Round to nearest 0.1 µL.

<sup>[2]</sup> Replace with Nuclease-free Water if performing endpoint PCR.

4. Remove the film seal from the sample plate, then add 30 µL Amplification Master Mix to each well (45-µL final volume).

5. Pierce the foil above the desired well of the Barcode Plate (see plate map) with a pipette tip. With a new tip, withdraw 5 µL of the Ion SingleSeq™ Barcode Adapter, then add to the appropriate sample (50-µL final volume). Repeat for each sample. Each Barcode Adapter is single-use only.

**Note:** To avoid sample misidentification later, be careful to track the correspondence of each sample and its Barcode Adapter.

**Note:** See “About the Ion SingleSeq™ Barcode Set 1” on page 12 for barcode plate handling guidelines.

	1	2	3	4	5	6	7	8	9	10	11	12
A	1	9	17									
B	2	10	18									
C	3	11	19									
D	4	12	20									
E	5	13	21									
F	6	14	22									
G	7	15	23									
H	8	16	24									

Barcode Plate containing Ion SingleSeq™ Barcode Set 1. Barcode Adapters are loaded in wells A1 through H3.

6. Adjust a pipettor to 30 µL, then mix the samples by pipetting up and down, using a new tip for each sample. Do not introduce bubbles into the samples.  
**Note:** If you used a non-optical reaction plate or PCR tube strip for cell lysis and pre-amplification, you can transfer samples to an optical plate at this point to perform real-time PCR amplification.
7. Apply a new film seal to the plate and briefly centrifuge to collect liquid at the bottom of the wells.
8. Cycle the samples in the thermal cycler using the following program:

Step	Temperature	Time	Number of cycles
1	95°C	3 minutes	1
2	95°C	20 seconds	4
	50°C	25 seconds	
	72°C	40 seconds	
3	95°C	20 seconds	12 <sup>[1]</sup>
	72°C	55 seconds	
4	4°C	Hold	1

<sup>[1]</sup> Acquire fluorescence data at this step if monitoring amplification in real-time.

9. Remove the plate, centrifuge at 1,000 × g for 30 seconds to collect liquid at the bottom of the wells, and place on ice or in a cold block.

Proceed to "Library pooling, purification, and quantification".

**STOPPING POINT** Sample libraries can be stored at –30°C to –10°C before proceeding to the next step.

## Library pooling, purification, and quantification

**Note:** We recommend that you perform library pooling, purification, and quantification in one session. Do not store library pools before quantification.

### Materials required

- 70% Ethanol - freshly prepared
- AMPure™ XP Reagent - warmed to room temperature
- DynaMag™-2 Magnet
- Low TE (10 mM Tris-HCl, 0.1 mM EDTA, pH 8.0)
- Qubit™ dsDNA HS Assay Kit
- Qubit™ 2.0 or Qubit™ 3.0 Fluorometer
- Qubit™ Assay Tubes

### Pool the libraries

Pool the libraries according to the PCR method you used to amplify the libraries. Before pooling, vortex the amplification reactions to mix, then centrifuge briefly to collect contents at the bottom of the wells or tubes.

If you used	Action
<b>Standard endpoint PCR amplification</b>	Add 5 µL of each library to a new 0.2-mL tube to create an equi-volume pool. Vortex the tube to mix and pulse-centrifuge to collect contents at the bottom of the tube.
<b>Real-time PCR amplification</b>	<p>Calculate the median <math>C_t</math> value of the libraries you want to pool and adjust the volumes of libraries deviating from the median <math>C_t</math> following these guidelines:</p> <ul style="list-style-type: none"> <li>• Libraries that fall within 1 <math>C_t</math> of the median should be added at the normal volume (5 µL).</li> <li>• Libraries that fall &gt;3 <math>C_t</math>s <b>later</b> than the median should not be included in the pool. It is likely that the input material was not present in the tube or well or was absorbed to the tube side.</li> <li>• Libraries that fall <math>\geq 1</math> but <math>\leq 3</math> <math>C_t</math>s <b>later</b> than the median should be added at 2X the normal volume (10 µL).</li> <li>• Do not adjust more than 25% of the libraries in a pool (excluding libraries &gt;3 <math>C_t</math>s later than median). If more than 25% of libraries fall 1 or more <math>C_t</math>s away from the median, adjust libraries with the most extreme <math>C_t</math> deviations until the 25% threshold is reached.</li> </ul>

**Note:** When pooling fewer than 8 libraries, the pool volume drops below 40 µL. Add Nuclease-free Water to bring the final volume to 40 µL before library pool purification.

## Purify the library pool

**Note:** Purify the *entire volume* of the library pool. For a pool of fewer than 8 libraries, bring the volume to 40 µL with Nuclease-free Water.

1. Heat the library pool in a thermal cycler using the following program:

Step	Temperature	Time	Number of cycles
1	70°C	2 minutes	1
2	22°C	Hold	1

2. Pulse-centrifuge the tube to collect contents, then transfer the heated library pool to a new 1.5-mL Eppendorf LoBind™ Tube.
3. Total the volumes of libraries and water added to the pool, then add an equal volume of AMPure™ XP Reagent beads.
4. Vortex briefly, pulse-centrifuge the tube to collect contents, then incubate for 5 minutes at room temperature.
5. Place the tube in the DynaMag™-2 magnet, then wait 5 minutes for beads to aggregate to the side of the tube.
6. Aspirate the supernatant carefully, then discard.
7. Wash beads with 150 µL of 70% ethanol while the tube is still on the magnet.
8. Incubate for 30 seconds.
9. Aspirate, then discard the wash solution.
10. Repeat steps 7–9, thoroughly removing all ethanol after the second wash.
11. Allow the beads to dry at room temperature for 3–4 minutes with the tube on the magnet.
12. Remove the tube from the magnet, then resuspend beads in 200 µL of Low TE by pipetting up and down.
13. Incubate the tube at room temperature for 1 minute.
14. Place the tube in the DynaMag™-2 magnet, then wait 2–3 minutes for beads to aggregate to the side of the tube.
15. Transfer 195 µL of the supernatant containing the purified library pool to a new 1.5-mL Eppendorf LoBind™ Tube and place on ice. Avoid carryover of beads.

## Quantify the library pool

Quantify the Ion SingleSeq™ library pool with the Qubit™ dsDNA HS (High Sensitivity) Assay Kit. See the *Qubit™ dsDNA HS Assay Kits User Guide* (Pub. No. MAN0002326) for more information.

1. Prepare a Qubit™ working solution by diluting the Qubit™ dsDNA HS Reagent (Component A) 1:200 in Qubit™ dsDNA HS Buffer (Component B) in a plastic tube.

Prepare sufficient working solution to quantify one or more purified pools plus two standards (volume required = (number of pools + 2) × 200 µL × 1.1).

---

**IMPORTANT!** Do not prepare the working solution in a glass container.

---

2. Prepare the standards:
  - a. Add 190 µL Qubit™ working solution to two labeled Qubit™ Assay Tubes used for standards.
  - b. Add 10 µL of each Qubit™ standard (Components C and D) to the appropriate tube.
  - c. Mix by vortexing 2–3 seconds. Do not create bubbles.
3. Prepare the unknown:
  - a. Aliquot 198 µL Qubit™ working solution to labeled Qubit™ Assay Tubes used for samples.
  - b. Add 2 µL of the library pool to the appropriate sample tube.
  - c. Mix by vortexing 2–3 seconds. Do not create bubbles.
4. Incubate the tubes in the dark at room temperature for 2 minutes.
5. Measure standards first to generate a standard curve, then measure the library pool. For the Qubit™ 2.0 instrument:
  - a. For the library pool: on the Qubit™ instrument, press **Calculate Stock Conc..**
  - b. Turn roller wheel to 2 (µLs original sample added to reagent).
  - c. Adjust units to ng/µL, then record this value.
6. Convert ng/µL to nM by multiplying the ng/µL values obtained in step 5 by 6.06 nmol/mg.

**Example:** Library pool concentration is 10 ng/µL

10 ng/µL (or 10 mg/L) × 6.061 nmol/mg = 60.6 nmol/L = 60.6 nM

**Note:** For DNA segments of 250-bp average length:

MW = (250 bp × 660,000 mg/mol/bp) = 1.65 × 10<sup>8</sup> mg/mol

Inverting 1.65 × 10<sup>8</sup> mg/mol and multiplying by 1 × 10<sup>9</sup> nmol/mol gives 6.06 nmol/mg.



7. Dilute the library pool to 1 nM.

In the preceding example, add 298  $\mu\text{L}$  of Low TE to 5  $\mu\text{L}$  of 60.6-nM pooled library stock.

**Note:** Diluted library pools and non-diluted library pool stock solutions can be stored for 1 week at 4°C. After 1 week, we recommend that you thaw the individual Ion SingleSeq™ libraries and repeat the pooling, purification, quantification, and dilution steps starting from “Pool the libraries” on page 30.



# Prepare template-positive Ion PGM<sup>™</sup> Template IA ISPs

■ Materials required .....	34
■ Isothermal amplification (IA) reaction procedural guidelines .....	35
■ Before you begin .....	35
■ Prepare the final library pool dilution .....	36
■ Perform the IA reaction .....	36
■ Recover the template-positive ISPs .....	38

## Materials required

### **Provided in the Ion PGM<sup>™</sup> Template IA Reagents 500 (Part No. A24619):**

- Ion PGM<sup>™</sup> Template IA Pellet (one pellet per reaction)
- Ion PGM<sup>™</sup> Template IA ISP Dilution Buffer
- Ion PGM<sup>™</sup> Template IA Start Solution

### **Provided in the Ion PGM<sup>™</sup> Template IA Reactions 500 (Part No. A24620):**

- Ion PGM<sup>™</sup> Template IA Ion Sphere<sup>™</sup> Particles (ISPs)
- Ion PGM<sup>™</sup> Template IA Primer Mix S
- Ion PGM<sup>™</sup> Template IA Rehydration Buffer

### **Provided in the Ion PGM<sup>™</sup> Template IA Solutions 500 (Part No. A24621):**

- Ion PGM<sup>™</sup> Template IA Stop Solution
- Ion PGM<sup>™</sup> Template IA Recovery Solution
- Ion PGM<sup>™</sup> Template IA Wash Solution

### **Other Materials and equipment**

- Nuclease-free Water
- 2-mL Eppendorf LoBind<sup>™</sup> Tubes
- Vortexer
- Microcentrifuge
- Low retention barrier pipette tips
- Pipettors
- Heat block set to 40°C (with water added to wells)

## Isothermal amplification (IA) reaction procedural guidelines

### Guidelines for preventing contamination

The IA reaction is highly sensitive to contaminating DNA. Follow the guidelines below to prevent introduction and carryover of contaminating DNA sequences in the work area.

- A primary source of contamination is spurious DNA fragments from previous sample processing steps. Do not introduce amplified DNA into the work area.
- To prevent carryover contamination, wipe down pipettors, hood and other work surfaces with a 70% ethanol-moistened Kimwipes™ disposable wipe before each experiment.
- After performing the reaction, rinse tube racks with Milli-Q™-quality water and dry.
- If possible, perform Ion PGM™ Template IA ISP preparation/reaction assembly and amplification/ISP recovery in two rooms: a pre-PCR clean room for ISP preparation/reaction assembly and a “dirty” post-PCR room for amplification and ISP recovery.
- If two rooms are used, ISP preparation and reaction assembly in a laminar flow hood is recommended but not required.
- If the entire procedure needs to be performed in one room, prepare the Ion PGM™ Template IA ISPs in a clean pre-PCR laminar flow hood.

### IA reaction guidelines

- Use Ion PGM™ Template IA Primer Mix S (black cap) in the Ion ReproSeq™ PGS procedure. Do not use Primer Mix L (blue cap).
- To achieve optimal results, do not agitate tubes after the pulse-centrifugation step that follows addition of the Ion PGM™ Template IA Start Solution to the IA reactions.
- The length of the IA reaction (25 minutes at 40°C) is important. Follow these guidelines if you are performing multiple reactions:
  - Limit the number of reactions you perform at the same time to 4.
  - If you need to perform more than 4 reactions, stagger the start and termination of each additional reaction by 5 minutes.

## Before you begin

1. Set a heat block to 40°C, with water added to wells.
2. Thaw the Ion PGM™ Template IA Primer Mix S (black cap), and keep it and the Ion PGM™ Template IA Start Solution (purple cap) on ice while setting up the reaction.
3. Thaw the Ion PGM™ Template IA ISPs (orange cap) and Ion PGM™ Template IA Rehydration Buffer (white cap) and keep on ice while setting up the reaction.

## Prepare the final library pool dilution

1. Dilute the 1-nM library pool prepared in the previous chapter 1:100 with Nuclease-free Water to generate a 10-pM library pool.
2. Transfer 50 µL of the 10-pM library pool to a new 0.2-mL PCR tube.
3. Heat the tube in a thermal cycler using the following program:

Step	Temperature	Time	Number of cycles
1	70°C	2 minutes	1
2	4°C	Hold	1

4. Remove the tube from the thermal cycler and place on ice.

## Perform the IA reaction

1. Prepare Templating Solution in a 2-mL Eppendorf LoBind™ Tube on ice (or a cold block):

Templating Solution

Order of addition	Component	Volume per reaction
1	Ion PGM™ Template IA ISP Dilution Buffer (yellow cap)	122 µL
2	Ion PGM™ Template IA Primer Mix S <sup>[1]</sup> (black cap)	8 µL
3	Ion PGM™ Template IA ISPs <sup>[2]</sup> (orange cap)	21 µL
4	Library pool (10 pM) <sup>[3]</sup>	18 µL
—	<b>Total</b>	<b>169 µL</b>

<sup>[1]</sup> Do not use Primer Mix L (blue cap).

<sup>[2]</sup> Vortex 30 seconds at maximum speed to resuspend immediately before addition.

<sup>[3]</sup> From step 4 of the previous section.

2. Vortex the tube containing the Templating Solution for 2 seconds at maximum setting to mix, pulse-centrifuge, then return the tube to ice.
3. Invert the Ion PGM™ Template IA Rehydration Buffer (white cap) 3 times to mix, then use 720 µL to rehydrate the Ion PGM™ Template IA Pellet. Vortex for 2 seconds at maximum setting, then pulse-centrifuge to collect contents at the bottom of the tube. Place the rehydrated pellet on ice or a cold block.

**Note:** The rehydrated Ion PGM™ Template IA Pellet is opaque.

4. Transfer the rehydrated Ion PGM™ Template IA Pellet to Templating Solution on ice, vortex for 2 seconds at maximum setting, then pulse-centrifuge.

---

**IMPORTANT!** The rehydrated pellet solution is viscous. Ensure that you transfer the entire volume by pulse-centrifuging the rehydration tube after transfer and pipetting any remaining volume into the Templating Solution.

---

5. Invert the Ion PGM™ Template IA Start Solution (purple cap) 3 times to mix, then add 300 µL to the Template/IA Solution using the reverse pipetting technique.

**Note:** If you are setting up more than one IA reaction, follow steps 5 through 7 for each reaction before starting the next reaction.

To use the reverse pipetting technique, perform the following steps:

- a. Set a 1-mL pipette to 300 µL.
  - b. Press the pipette knob to the second stop, then dip the tip into the Start Solution.
  - c. Slowly release the pipette knob until it returns to the starting position. Allow 10 seconds for Start Solution to be fully drawn up into the tip (Fig. 5a).
  - d. Dispense the solution into the Templating Solution tube by gently pressing the pipette knob to the first stop point only. Wait at least 5 seconds until the liquid in the pipette tip stops moving. Some liquid remains in the tip (Fig. 5b).
  - e. Withdraw the tip from the tube. If any liquid adheres to the outer surface of the tip, touch the tip to the inner wall of the tube to transfer the liquid to the tube.
6. Vortex the tube 10 times in 1-second pulses at the maximum vortexer setting. Invert the tube, then repeat the ten 1-second pulses.
  7. Pulse-centrifuge the tube to collect contents, then immediately place the tube on ice.




---

**IMPORTANT!** Handle the tube gently after centrifuging. To achieve optimum results, do not agitate the tubes from this point on.

---

8. Start the IA reaction by gently placing the tube in the 40°C heat block. Ensure that the tube is immersed in water.
9. Incubate the IA reaction for 25 minutes at 40°C.

## Recover the template-positive ISPs

1. Stop the IA reaction by removing the tube from the heat block and adding 650 µL of Ion PGM™ Template IA Stop Solution.
2. Vortex the tube well to mix contents thoroughly, then centrifuge the tube at  $7,500 \times g$  for 3 minutes.
3. Aspirate, then discard the supernatant, being careful not to disturb the pellet. Leave ~100 µL in the tube.
4. Resuspend the pellet in 1 mL Ion PGM™ Template IA Recovery Solution.
  - a. Pipette up and down to resuspend the pellet.
  - b. Add an additional 700 µL Ion PGM™ Template IA Recovery Solution, then vortex thoroughly.
5. Incubate for 5 minutes with vortexing 5 seconds every minute.
6. Centrifuge for 3 minutes at  $12,000 \times g$ .
7. Immediately remove, then discard all of the supernatant without disturbing the ISP pellet. Remove any bubbles before removing the bulk of the liquid to avoid frothing in subsequent steps.

**Note:** The ISPs form a glassy pellet that is barely visible. Note the orientation of the tube in the centrifuge so that the position of the ISP pellet is known. The supernatant must be removed immediately to minimize the resuspension of ISPs.
8. Add 100 µL of the Ion PGM™ Template IA Wash Solution to the ISP pellet.
9. Resuspend the templated ISPs completely by vortexing for 4 seconds at maximum speed, then pipet the ISP suspension up and down 4 times. Proceed to Chapter 5, “Enrich the template-positive Ion PGM™ Template IA ISPs”.

---

**STOPPING POINT** Store templated ISPs in Ion PGM™ Template IA Wash Solution at 4°C for up to one week.

---



# Enrich the template-positive Ion PGM™ Template IA ISPs

■ Materials required .....	39
■ Determine if a residual volume test is necessary .....	40
■ Prepare reagents then fill the 8-well strip .....	41
■ Prepare the Ion OneTouch™ ES .....	44
■ Perform the Ion OneTouch™ ES run .....	46
■ Sequence or store the template-positive ISPs .....	48
■ Perform quality control on enriched ISPs .....	48

## Materials required

### Provided in Ion PGM™ Template IA Solutions 500 (Part No. A24621):

- Ion OneTouch™ Template IA Wash Solution
- MyOne™ Beads Wash Solution
- Tween™ Solution
- Neutralization Solution

### Provided in Ion PGM™ Template IA Supplies 500 (Part No. A24618):

- 8-well strip
- Eppendorf™ LoRetention Dualfilter Tips (P300)

### Other Materials and Equipment:

- Ion OneTouch™ ES Instrument
- Ion PGM™ Enrichment Beads (Cat. No. 4478525; Dynabeads™ MyOne™ Streptavidin C1 Beads)
- 1.5-mL Eppendorf LoBind™ Tubes
- 0.2-mL PCR tubes
- Nuclease-free Water
- 1 M NaOH
- Pipettes
- Vortexer
- DynaMag™-2 magnet
- Microcentrifuge

## Determine if a residual volume test is necessary

**IMPORTANT!** Ensure that the AC line voltage module is installed correctly into the Ion OneTouch™ ES Instrument. See the *Ion OneTouch™ 2 System User Guide* (Pub. No. MAN0014388) for information regarding instrument setup, calibration, and maintenance.

To determine if a residual volume test is necessary, follow these guidelines:

Condition	Action
First use of the instrument and during monthly maintenance	Perform a residual volume test (see "Ion OneTouch™ ES Instrument installation, setup, and maintenance" in the <i>Ion OneTouch™ 2 System User Guide</i> Pub. No. MAN0014388).
Routine use and residual volume in Well 1 and Well 8 is >5.0 µL	
Routine use and residual volume in Well 1 and Well 8 is ≤5.0 µL	Operate the instrument without performing the residual volume test. Proceed to "Prepare reagents then fill the 8-well strip" on page 41.



## Prepare reagents then fill the 8-well strip

### Prepare Melt-Off Solution

Prepare fresh Melt-Off Solution by combining the components in the following order:

Order	Component	Volume
1	Tween™ Solution	280 µL
2	1 M NaOH	40 µL
—	<b>Total</b>	<b>320 µL</b>

**IMPORTANT!** Prepare Melt-Off Solution as needed, but appropriately dispose of the solution after 1 day.

The final composition of the Melt-Off Solution is 125 mM NaOH and 0.1% Tween™ 20 detergent.

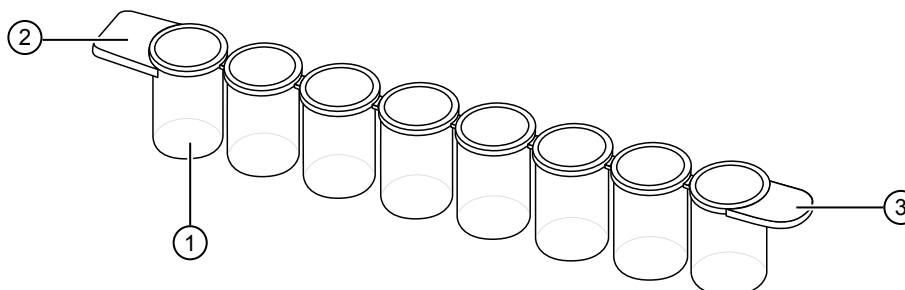
### Wash and resuspend the Dynabeads™ MyOne™ Streptavidin C1 Beads

1. Vortex the tube of Dynabeads™ MyOne™ Streptavidin C1 Beads for 30 seconds to resuspend the beads thoroughly, then centrifuge the tube for 2 seconds.
2. Open the tube, then use a new tip to pipet the dark pellet of beads up and down until the pellet disperses. Immediately proceed to the next step.
3. Transfer 13 µL of Dynabeads™ MyOne™ Streptavidin C1 Beads to a new 1.5-mL Eppendorf LoBind™ Tube.
4. Place the tube on a magnet such as a DynaMag™-2 magnet for 2 minutes, then carefully remove and discard the supernatant without disturbing the pellet of Dynabeads™ MyOne™ Streptavidin C1 Beads.
5. Add 130 µL of MyOne™ Beads Wash Solution to the Dynabeads™ MyOne™ Streptavidin C1 Beads.
6. Remove the tube from the magnet, vortex the tube for 30 seconds, then centrifuge for 2 seconds.

## Fill the 8-well strip

**Note:** If you stored template-positive ISPs at 2°C to 8°C, vortex the tube to resuspend the ISPs and pulse-centrifuge to collect contents. Pipet the solution up and down to resuspend the Ion PGM™ Template IA ISPs and transfer to Well 1 of the 8-well strip.

1. Add the entire volume (~100 µL) of template-positive ISPs from the amplification reaction into Well 1 of the 8-well strip. Well 1 with the ISPs is on the left:



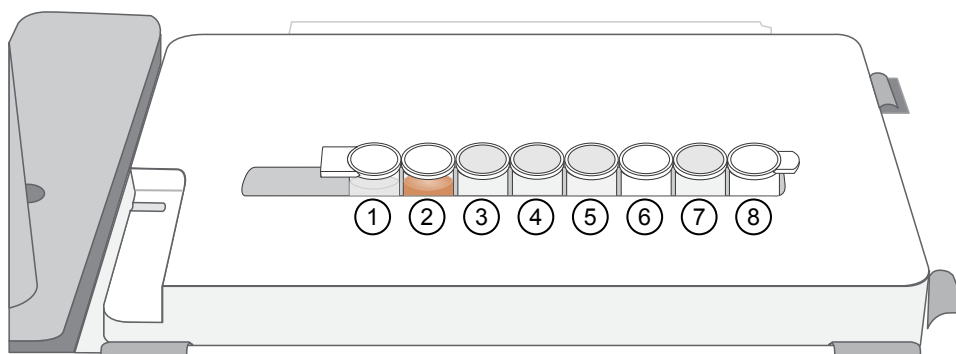
- ① Well 1
- ② Square-shaped tab
- ③ Rounded tab

2. If you have not already assessed the quality of the unenriched, template-positive ISPs, use the following method:

Quality assessment by	Action
Guava™ easyCyte™ 5 Flow Cytometer	Transfer a 1.0-µL aliquot of the unenriched ISPs to a 1.5-mL Eppendorf LoBind™ Tube. See the <i>Ion Sphere™ Particles (ISPs) Quality Assessment Using the Guava™ easyCyte™ 5 Flow Cytometer User Bulletin</i> (Pub. No. MAN0015799), available at: <a href="http://tools.thermofisher.com/content/sfs/manuals/MAN0015799.pdf">http://tools.thermofisher.com/content/sfs/manuals/MAN0015799.pdf</a>
Demonstrated protocol: Quality assessment by the Applied Biosystems™ Attune™ Acoustic Focusing Cytometer	Transfer a 1.0-µL aliquot of the unenriched ISPs to a 1.5-mL microcentrifuge tube. Put the sample on ice, then see the <i>Demonstrated Protocol: Ion Sphere™ Particles (ISPs) Quality Assessment using the Applied Biosystems™ Attune™ Acoustic Focusing Cytometer User Bulletin</i> (Pub. No. 4477181), available at: <a href="http://tools.thermofisher.com/content/sfs/manuals/4477181A.pdf">http://tools.thermofisher.com/content/sfs/manuals/4477181A.pdf</a>

3. Fill the remaining wells in the 8-well strip as follows (see the following figure):

Well number	Reagent to dispense in well
Well 1 (well nearest to the square-shaped tab)	Entire template-positive ISP sample (100 µL; prepared in step 1 of this procedure)
Well 2	130 µL of Dynabeads™ MyOne™ Streptavidin C1 Beads resuspended in MyOne™ Beads Wash Solution [prepared in “Wash and resuspend the Dynabeads™ MyOne™ Streptavidin C1 Beads” on page 41]
Well 3	300 µL of Ion PGM™ Template IA Wash Solution
Well 4	300 µL of Ion PGM™ Template IA Wash Solution
Well 5	300 µL of Ion PGM™ Template IA Wash Solution
Well 6	Empty
Well 7	300 µL of freshly-prepared Melt-Off Solution (prepared in “Prepare Melt-Off Solution” on page 41)
Well 8	Empty

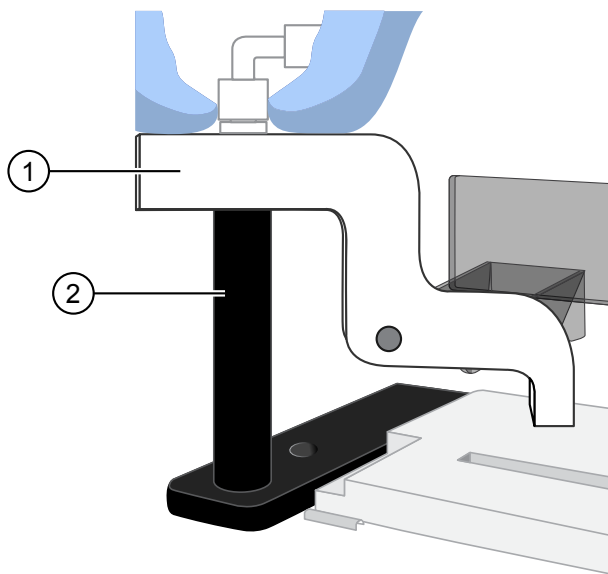


4. Ensure that the square-shaped tab is on the left, then insert the filled 8-well strip with the 8-well strip pushed all the way to the right end of the slot of the Tray.

## Prepare the Ion OneTouch™ ES

Before every enrichment performed on the Ion OneTouch™ ES Instrument, install a new Eppendorf™ LoRetention Dualfilter P300 pipette tip.

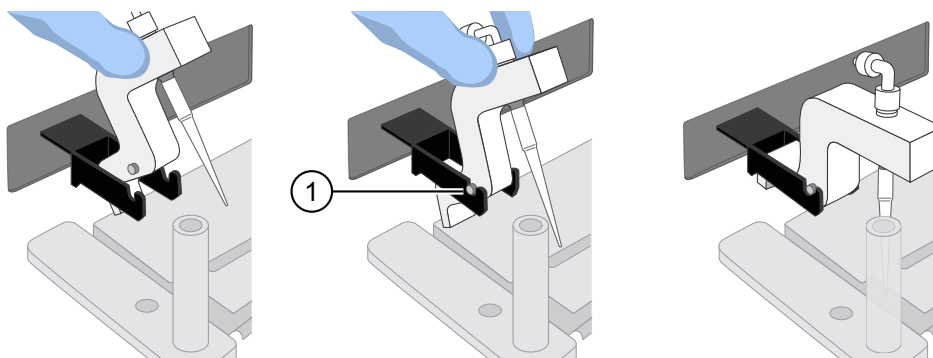
1. Place a new tip in the Tip Loader. Remove the Tip Arm from the cradle, then align the metal fitting of the Tip Arm with the tip.
2. Keeping the fitting on the Tip Arm vertical, firmly press the Tip Arm down onto the new tip until the Tip Arm meets the Tip Loader. Hold the Tip Arm to the Tip Loader for ~1 second to ensure proper installation of the tip.



- ① Tip Arm  
② Tip Loader

3. Lift the Tip Arm *straight* up to pull the installed tip from the Tip Loader tube.
4. Return the Tip Arm to the cradle.
  - a. Tilt the Tip Arm back (below left), then align the pins with the round notches in the cradle (below center).
  - b. Lower the Tip Arm into position (below center).

- c. Move the Tip Arm forward into the working position (below right).



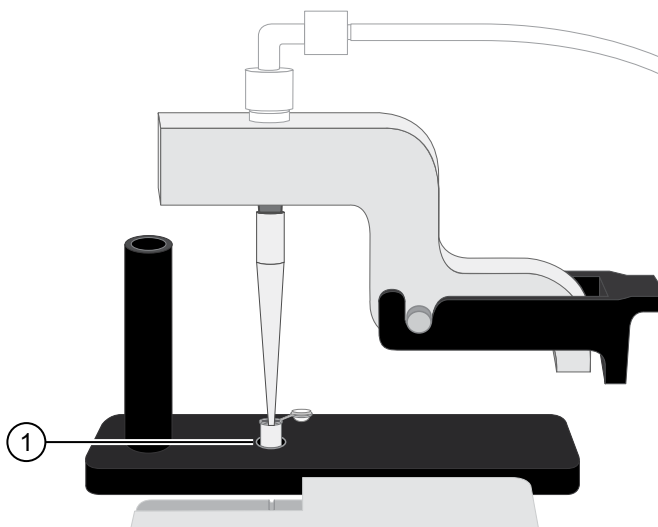
- ① Tip Arm pins resting in the notches in the cradle

---

**IMPORTANT!** Ensure that the back/bottom end of the Tip Arm is not resting on top of the thumb screw, causing the Tip Arm to tilt forward.

---

5. Add 10  $\mu$ L of Neutralization Solution to a new 0.2-mL PCR tube.
6. Insert the open 0.2-mL PCR tube containing Neutralization Solution into the hole in the base of the Tip Loader, as shown in the following figure.



- ① 0.2-mL PCR tube placed in hole at base of Tip Loader

## Perform the Ion OneTouch™ ES run

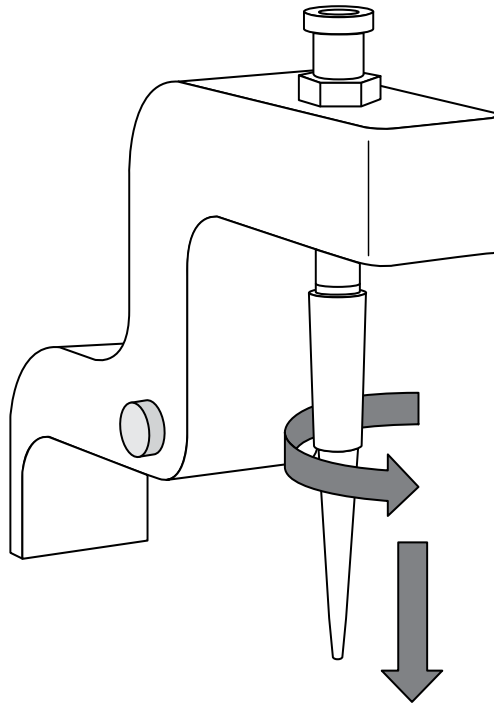
Ensure that a new tip and opened 0.2-mL PCR tube with the Neutralization Solution have been loaded. Ensure that Well 1 (ISP sample) is the left-most well and that the 8-well strip is pushed to the far-right position in the slot.

1. Pipet the contents of Well 2 up and down to resuspend the beads before starting the run. Do not introduce bubbles into the solution.
2. If needed, power on the Ion OneTouch™ ES, then wait for the instrument to initialize. The screen displays “rdy”. The Tip Arm performs a series of initialization movements and returns to the home position (~5 seconds).
3. Press **Start/Stop**. The screen displays “run” during the run. The run takes ~35 minutes.

**Note:** If necessary to stop a run, press **Start/Stop**. The instrument completes the current step, then stops the run and displays “End”. Press **Start/Stop** again to return the Tip Arm to the home position. It is not possible to restart (where you left off) after stopping a run.

4. At the end of the run, the instrument displays “End” and beeps every 60 seconds. Press the **Start/Stop** button to silence this alarm, then reset the Ion OneTouch™ ES for the next run. The instrument can be left on between runs.
5. *Immediately after the run*, securely close and remove the PCR tube containing the enriched ISPs.
6. Mix the contents of the PCR tube by gently inverting the tube five times.  
Ensure that the 0.2-mL PCR tube has >200 µL of solution containing the enriched ISPs. After a successful run on the instrument, the sample is in ~230 µL of Melt-Off Solution, Ion PGM™ Template IA Wash Solution, and Neutralization Solution. If the tube has <<200 µL of solution containing the enriched ISPs, contact Technical Support.

7. Remove the used tip: While you are standing above the Tip Arm, and with the Tip Arm in its cradle, twist the tip counterclockwise, then pull it downward to remove. Discard the tip:



---

**IMPORTANT!** Improper removal of tips can loosen the metal tip adapter fitting on the Tip Arm and affect instrument operation.

---

8. Remove, then discard the used 8-well strip.

## Sequence or store the template-positive ISPs

- Sequence using the Ion PGM™ Hi-Q™ Sequencing Kit, bundled in the Ion ReproSeq™ PGS Kit. Proceed to Chapter 6, “Create a Planned Run”.
- or
- Store the enriched ISPs at 2°C to 8°C for up to 3 days.

## Perform quality control on enriched ISPs

You can determine the enrichment efficiency using one of the following methods:

Quality assessment by	Action
Guava™ easyCyte™ 5 Flow Cytometer	Transfer a 1.0-μL aliquot of the enriched ISPs to a 1.5-mL Eppendorf LoBind™ Tube. Refer to the <i>Ion Sphere™ Particles (ISPs) Quality Assessment Using the Guava™ easyCyte™ 5 Flow Cytometer User Bulletin</i> (Pub. No. MAN0015799), available at: <a href="http://tools.thermofisher.com/content/sfs/manuals/MAN0015799.pdf">http://tools.thermofisher.com/content/sfs/manuals/MAN0015799.pdf</a>
Attune™ NxT Acoustic Focusing Cytometer	Transfer a 1.0-μL aliquot of the enriched ISPs to a 1.5-mL Eppendorf LoBind™ Tube. Put the sample on ice, then refer to <i>Demonstrated Protocol: Ion Sphere™ Particles (ISPs) Quality Assessment using the Applied Biosystems™ Attune™ Acoustic Focusing Cytometer User Bulletin</i> (Pub. No. 4477181), available at: <a href="http://tools.thermofisher.com/content/sfs/manuals/4477181A.pdf">http://tools.thermofisher.com/content/sfs/manuals/4477181A.pdf</a>





# Create a Planned Run

■ About Planned Runs .....	49
■ Create a Planned Run .....	50
■ Planned Run wizard: key fields .....	53

---

**IMPORTANT!** This sequencing kit is compatible with Torrent Suite™ Software v5.0 and later. Before proceeding, check for updates to the Torrent Suite™ and sequencing instrument software, and install the updates if available.

---

## About Planned Runs

Planned Runs contain all the settings used in a sequencing run, including number of flows, kit types, barcodes used (if any), run type (e.g., DNA, RNA, amplicons), and reference file (if any). They provide a fast and convenient way to set up and organize your runs.

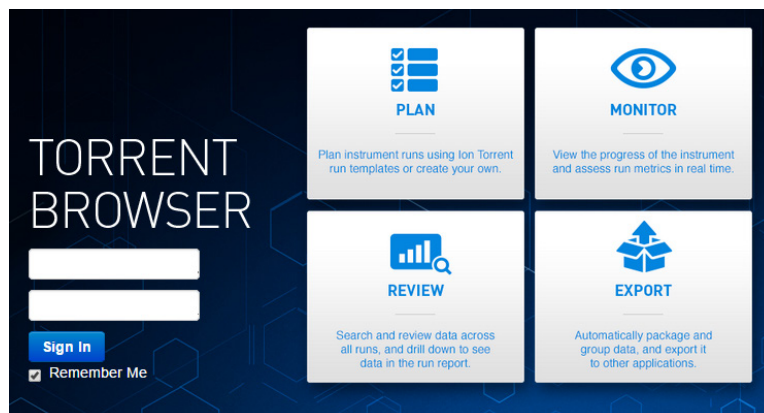
You create a Planned Run using Torrent Browser on the Torrent Server connected to your sequencer, and then select the appropriate plan in the **Select Planned Run** screen of the sequencer touchscreen when you start the run.

You can also create a Planned Run on one Torrent Server and then transfer it to another server for sequencing. See the *Ion PGM™ Hi-Q™ Sequencing User Guide*, (Pub. No. MAN0009816; Appendix B), for more information.

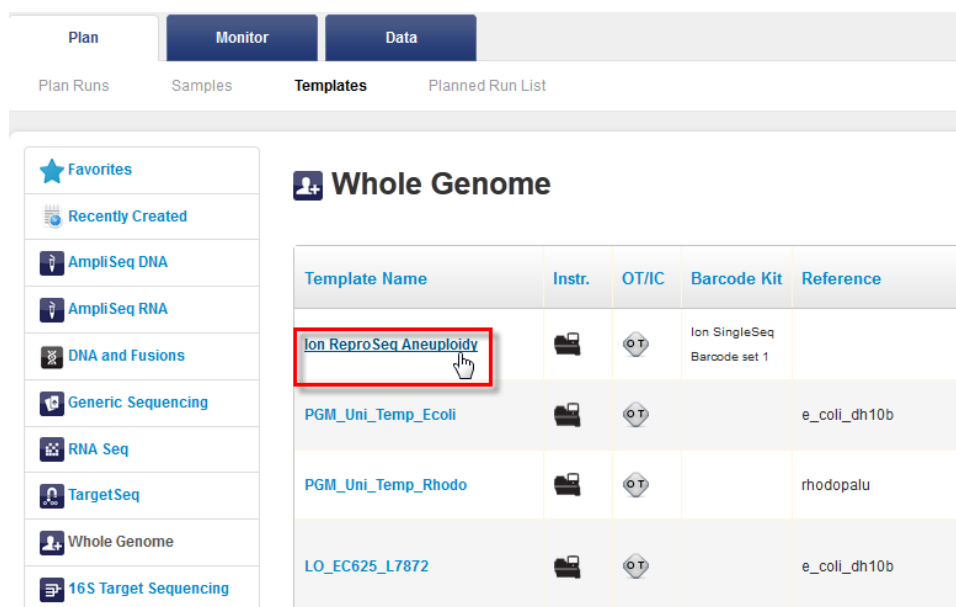
**Note:** For additional information, see the *Torrent Suite™ Software Help Guide* (Pub. No. MAN0014688), available at: <https://tools.thermofisher.com/content/sfs/manuals/MAN0014668.pdf>.

## Create a Planned Run

1. Open the Torrent Browser on the Torrent Server connected to your sequencer.



2. Select the **Plan** tab, select **Templates**, then click **Whole Genome** from the list of applications on the left side of the page.
3. Select **Ion ReproSeq Aneuploidy** from the list of Planned Run templates.



4. The Ion ReproSeq™ Aneuploidy template will auto-populate settings in the Application, Kits, and Plugins tabs in the Planned Run wizard appropriately. Key fields are described in “Planned Run wizard: key fields” on page 53.

**Note:** Select the appropriate chip type you are using in the **Kits** tab. Chip type is set to Ion 318™ Chip v2 by default.

In the **Plan** tab,

- a. Enter a new Run Plan Name, if desired.
- b. Select **hg19(Homo sapiens)** from the Reference Library dropdown list.

- c. Select **None** from the Target Regions and Hotspot Regions dropdown lists.
- d. Enter the number of barcodes you are using in your Sample Set.
- e. Select a Barcode from the dropdown list for each sample, if different from the Barcode shown. Enter a unique, descriptive name, such as one including one or more sample-specific or grouping attributes you wish to track, for each Ion SingleSeq™ Barcode used. Avoid use of the default "Sample 1", "Sample 2", etc.

Template Name :  
Ion ReproSeq Aneuploidy

Run Plan Name (required) :  
Ion ReproSeq Aneuploidy

Analysis Parameters: ☒ Default (Recommended) ☐ Custom [Details +](#)

Default Reference & BED Files

Reference Library : hg19(Homo sapiens)

Target Regions: None

Hotspot Regions: None

☒ Use same reference & BED files for all barcodes

Number of barcodes : 24 [Save Samples Table](#) [Load Samples Table](#)

Sample Tube Label :

Chip ID :

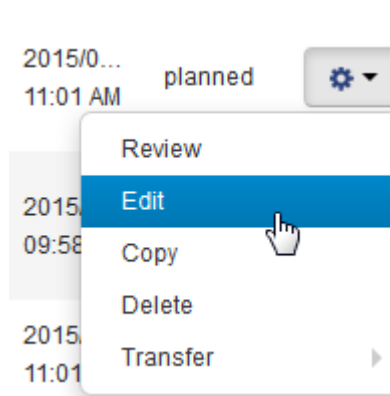
Enter a sample name for each barcode used (require at least one sample) [↓](#) [🗑](#) [👁](#) :

#	Barcode	Sample Name (required)	Sample ID	Sample Description
1	SingleSeq_001 (TAGGTGGTTC)	Sample 1		
2	SingleSeq_002 (TCTATTCGTC)	Sample 2		
3	SingleSeq_003 (TCGCAATTAC)	Sample 3		

**IMPORTANT!** We strongly recommend that you assign unique sample names for each barcode and experiment. This will enable easy identification and retrieval of results in Ion Reporter™ analyses based on experiment name, plate code, etc. Do not reuse sample names from experiment to experiment.

- f. Click **Plan Run** in the lower right corner to save the Planned Run. The run is listed on the Planned Runs page under the name you entered.

5. To set up auto-analysis using a custom Ion Reporter™ workflow, navigate to your Planned Run in the Planned Run list, click the Settings symbol ⚙️ to the right of your run, then select **Edit** from the dropdown list.



6. Click the **IonReporter** tab, then
  - a. Select your account.
  - b. Select the **ReproSeq Low-pass whole-genome aneuploidy** workflow.
7. Complete your selections in the Projects tab, then click **Next**.
8. Click **Update Plan** in the lower right corner of the page to save your edits.

## Planned Run wizard: key fields

Field name	Description
IonReporter	Select the account and select ReproSeq Low-pass whole-genome aneuploidy workflow from the Existing Workflow menu. To create a new workflow, click <b>Create New Workflow</b> .
Application	Select the sequencing application you are performing: <b>DNA Whole Genome</b>
Library Kit Type	Select <b>Ion SingleSeq Kit</b> .
Template Kit	Select <b>Ion PGM Template IA 500 Kit</b> .
Sequencing Kit	Select <b>Ion PGM™ Hi-Q™ Sequencing Kit</b> .
Flows	Enter the appropriate number of flows for the sequencing kit and read length: <b>250 flows</b>
Chip Type	Select the Ion chip type you are using.
Forward 3' Adapter	Select <b>Ion P1</b> .
Barcode Set	Select <b>Ion SingleSeq Barcode set 1</b> .
Project	Select or add a project within which to group your run data: Ion ReproSeq™ PGS
Run Plan Name	Enter a name for the Planned Run.
Reference Library	Select a reference library uploaded to the Torrent Server: <b>hg19</b>
Target Regions and Hotspot Regions	Set to "None" (default).
Enter a sample name	Enter a Sample Name, set relation to "Self", and assign unique Analysis IDs for each sample in the run (number of samples will change based on the number of barcodes selected). Avoid using the default names "Sample 1", "Sample 2", etc.
Monitoring Thresholds	Set thresholds for Bead Loading, Key Signal, and Usable Sequence. In the Torrent Browser <b>Monitor ▶ Runs in Progress</b> tab, an alert is displayed if the values for a run fall below the selected thresholds.



# Clean and initialize the Ion PGM™ Sequencer

■ Condition the Wash 2 Bottle for first use .....	54
■ Cleaning .....	54
■ Initialize the Ion PGM™ System .....	57

## Condition the Wash 2 Bottle for first use

New Wash 2 Bottles must be conditioned with Wash 2 Bottle Conditioning Solution for at least 8 hours before first use.

**Note:** If necessary, you can reuse an existing Wash 2 Bottle while you condition a new bottle. Bottles can be used for sequencing up to 40 times before they must be replaced.

To condition the Wash 2 Bottle:

1. Fill the bottle to the mold line with 18 MΩ water, add the entire container of Wash 2 Bottle Conditioning Solution, then cap the bottle and invert it 5 times to mix.
2. Allow the bottle to sit at room temperature for at least 8 hours and preferably overnight, then dispose of the contents. The bottle is now ready for use.

## Cleaning

### Materials required

- 18 MΩ water (e.g., the ELGA™ PURELAB™ Flex Water Purification System)
- Cleaning bottles and collection trays (provided with the Ion PGM™ System)
- Old chip that has been used for sequencing, marked for cleaning
- Used sipper tubes (from the previous run)
- Squirter bottle
- **Chlorite cleaning:** Ion PGM™ Cleaning Tablet (provided in the kit)
- **Chlorite cleaning:** 1 M NaOH, diluted fresh each week from 10 M NaOH
- **Chlorite cleaning:** Glass bottle (1 L)
- **Chlorite cleaning:** 0.22-μm or 0.45-μm vacuum filtration system and filters

## Cleaning schedule

The Ion PGM™ Sequencer requires cleaning with either 18-MΩ water or a chlorite solution every time the instrument is initialized.

Clean with	Schedule
18 MΩ water	<ul style="list-style-type: none"><li>• Daily, when instrument is in use (e.g., not necessary on weekends)</li><li>• After one or more runs totaling ≤1,100 flows</li><li>• If more than 27 hours but less than 48 hours have elapsed between the last cleaning/initialization and the start of a run</li><li>• If you cleaned with chlorite a week ago and have not used the instrument since then</li></ul>
Chlorite solution	<ul style="list-style-type: none"><li>• Once a week, unless the instrument has not been used since the last chlorite cleaning (in which case, clean with 18 MΩ water before using)</li><li>• If the instrument has been left with reagents for more than 48 hours (for example, over the weekend)</li></ul>

## Cleaning setup

**IMPORTANT!** For all the following steps, use 18 MΩ water directly from the purification system. Do not use water that has been collected or stored in any other containers.

- Remove any wash and reagent bottles that are attached to the Ion PGM™ System before cleaning.
- Do not remove old sippers before cleaning. The sippers are used as part of the cleaning procedure.
- Old chips that have been used for sequencing can be marked and used in the cleaning procedure.
- Wash bottles (250 mL and 2 L) provided as part of instrument installation can be marked and used for cleaning. After you have used the wash bottles provided with the sequencing kit for the specified number of runs, you can use them as extra cleaning bottles. Mark them for cleaning use only.

## 18 MΩ water cleaning

1. Empty any remaining solution from each cleaning bottle (two 250-mL bottles and one 2-L bottle) and rinse each bottle twice with ~100 mL of 18 MΩ water.
2. Press **Clean** on the touchscreen, and select the **18-MOhm water cleaning** checkbox. Press **Next**.
3. Using ungloved hands, secure a used chip designated for cleaning in the chip clamp.

**IMPORTANT!** Always make sure that both red rubber gasket port fittings are securely in place when securing chips with the chip clamp. Failure to do so can result in a spill hazard and instrument damage.

4. Remove all wash and reagent bottles attached to the instrument. Keep the sippers in place at all positions. Press **Next**.
5. Add 250 mL of 18 MΩ water to an empty 250-mL cleaning bottle.

6. Rinse the outside of the sipper tube in the W1 position on the instrument with a squirt bottle containing 18 MΩ water.
7. Attach the 250-mL bottle containing 18 MΩ water to the W1 position, ensuring that the W1 cap is screwed on tightly. Press **Next**.
8. Place the empty 2-L cleaning bottle in the W2 position and the empty 250-mL bottle in the W3 position, and insert the sippers into the bottles. Do not screw on the caps.
9. Place collection trays below the reagent sippers in the dNTP positions. Press **Next** to begin cleaning.
10. When cleaning is complete, remove the bottles and sippers from the W1, W2 and W3 positions. Leave the reagent sippers and collection trays in place. Press **Next** to return to the main menu and proceed to initialization.

## Chlorite cleaning

**Note:** Prepare a stock of 1 M NaOH each week by diluting 10 M NaOH with 18 MΩ water.

1. Empty any remaining solution from each cleaning bottle (two 250-mL bottles and one 2-L bottle), then rinse each bottle twice with ~100 mL of 18 MΩ water.
2. Fill a glass bottle with 1 L of 18 MΩ water, then add an Ion Cleaning tablet (chlorite tablet). Allow the tablet to dissolve completely (~10 minutes).
3. When the tablet has dissolved, add 1 mL of 1 M NaOH and filter the solution using a 0.22-μm or 0.45-μm filter. Use the chlorite solution within 2–3 hours. Discard any unused solution after this time.
4. Press **Clean** on the touchscreen, then select the **Chlorite cleaning** checkbox. Press **Next**.
5. Using ungloved hands, secure a used chip designated for cleaning in the chip clamp.

---

**IMPORTANT!** Always ensure that both red rubber gasket port fittings are securely in place when securing chips with the chip clamp. Failure to do so can result in a spill hazard and instrument damage.

---

6. Remove all wash and reagent bottles that are attached to the instrument. Keep the sippers in place at all positions. Press **Next**.
7. Add 250 mL of the filtered chlorite solution to an empty 250-mL cleaning bottle.
8. Rinse the outside of the sipper tube in the W1 position on the instrument with a squirt bottle containing 18 MΩ water.
9. Attach the 250-mL bottle with the filtered chlorite solution to the W1 position. Ensure that the W1 cap is tight. Press **Next**.
10. Place the empty 2-L cleaning bottle in the W2 position and the empty 250-mL bottle in the W3 position, then insert the sippers into the bottles. Do not screw on the caps.



11. Place collection trays below the reagent sippers in the dNTP positions. Press **Next** to start cleaning.
12. When prompted, remove the bottle containing the chlorite solution from the W1 position.
13. Rinse the outside of the W1 sipper tube with a squirt bottle containing 18 MΩ water.
14. Fill a clean 250-mL bottle with 250 mL of 18 MΩ water, then attach the bottle in the W1 position. Ensure the cap is tight. Press **Next** to start the water rinse.
15. When cleaning is complete, remove the bottles and sippers from the W1, W2 and W3 positions. Leave the reagent sippers and collection trays in place. Press **Next** to return to the main menu, then proceed to initialization.

## Initialize the Ion PGM™ System

Initialization takes ~1 hour. As part of the initialization process, first prepare the Wash and Reagent Bottles as described in this section.

### Materials required

#### Materials provided in the kit

- Ion PGM™ Hi-Q™ Sequencing dGTP
- Ion PGM™ Hi-Q™ Sequencing dCTP
- Ion PGM™ Hi-Q™ Sequencing dATP
- Ion PGM™ Hi-Q™ Sequencing dTTP
- Ion PGM™ Hi-Q™ Sequencing W2 Solution (stored protected from light)
- Ion PGM™ Sequencing W3 Solution
- Wash 1 and Wash 3 Bottles and sipper tubes
- Wash 2 Bottle and sipper tubes (bottle must be conditioned prior to first use, as described in “Condition the Wash 2 Bottle for first use” on page 54)
- Wash 2 Bottle Conditioning Solution
- Reagent Bottles and sipper tubes

#### Other materials and equipment

- Used chip (leave chip on the instrument during initialization)
- 18 MΩ water
- 100 mM NaOH (prepared daily)
- Ice
- 5-mL and 25-mL pipettes
- Filtered and unfiltered pipette tips and pipettes
- Vortex mixer
- Microcentrifuge
- **Optional:** Ion PGM™ Sequencing Sippers Kit (Cat. No. 4478682)

## Initialization guidelines

---

**IMPORTANT!** Handle nucleotides carefully to avoid cross-contamination. Always change gloves after removing used sipper tubes from the Ion PGM™ System to avoid cross contamination of the nucleotides. Also change gloves after handling concentrated dNTP stocks.

---

For each initialization, the first run should be started within 1 hour after initialization, and the last run must be started within 24 hours after initialization.

### Bottle usage

- Wash 2 Bottles can be used for up to 40 initializations, after which you can use them in the cleaning procedure.
- Wash 1 and Wash 3 Bottles can be used for up to 4 initializations, after which you can reuse them in the cleaning procedure.
- Replace the Reagent Bottles and sipper tubes every time you initialize.

## Before initialization

1. Remove the dNTP stock solutions from the freezer and begin thawing on ice.
2. Check the tank pressure for the nitrogen gas. When the tank pressure drops below 500 psi, change the tank.

## Prepare the Wash 2 Bottle

### Note:

- Do not remove the old sippers from the dNTP ports until instructed to do so.
- Load the bottles as quickly as possible to prevent atmospheric CO<sub>2</sub> from reducing the pH of the Wash 2 solution.
- For all the following steps, pour the 18 MΩ water directly from the purification system into the Wash 2 Bottle. Do not use water that has been collected or measured in any other containers.

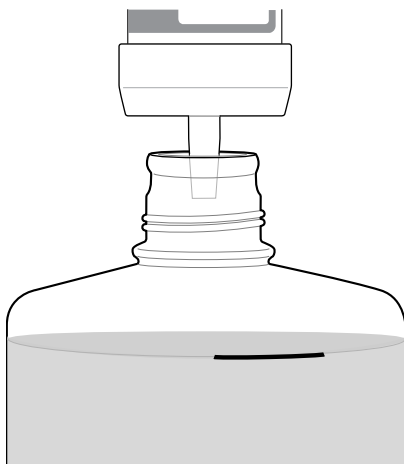
---

**IMPORTANT!** Do not let the new sippers touch any surfaces.

---

1. Rinse the Wash 2 Bottle (2 L) 3 times with 200 mL of 18 MΩ water.
2. Prepare 500 μL of 100 mM NaOH by diluting 50 μL of 1 M NaOH in 450 μL of nuclease-free water.

3. If your 18 MΩ water system has a spigot, extend it into **but not below** the neck of the Wash 2 Bottle. Otherwise, position the nozzle as close to the mouth of the bottle as possible.



**Note:** If your water system has a digital display, ensure it reads "18 MΩ" throughout filling the bottle. If not, see Appendix A, "Troubleshooting".

4. Fill the bottle to the mold line with 18 MΩ water. The volume of water is ~2 liters. (You can mark the mold line on the bottle for clarity.)

**Note:** If you are preparing bottles for multiple sequencers, cap each bottle immediately after filling, and leave capped until you are ready to add Ion PGM™ Hi-Q™ Sequencing W2 Solution.

5. Add the entire bottle of Ion PGM™ Hi-Q™ Sequencing W2 Solution to the Wash 2 Bottle.



**Note:** Keep the Ion PGM™ Hi-Q™ Sequencing W2 Solution bottle to scan the barcode during the initialization procedure.

- Using a P200 pipette, add 70 µL of 100 mM NaOH to the Wash 2 Bottle.

**Note:** Different sites can require adding different volumes of 100 mM NaOH. Some sites, for example, can require doubling the volume to 140 µL. See “Error message: Added too much W1 to W2” on page 98 for information on determining the volume of 100 mM NaOH to add.

- Cap the bottle and invert 5 times to mix, and immediately proceed through the remainder of the initialization procedure.

---

**IMPORTANT!** Do not store the mixed Wash 2 Bottle.

---

## Prepare the Wash 1 and Wash 3 Bottles

**Note:** For the following steps, label the Wash 1 and Wash 3 Bottles to avoid confusion.

- Rinse the Wash 1 and Wash 3 Bottles 3 times with 50 mL of 18 MΩ water.
- Wash 1 Bottle:** Add 350 µL of freshly prepared 100 mM NaOH to the Wash 1 Bottle, then cap the bottle.
- Wash 3 Bottle:** Add Ion PGM™ Sequencing W3 Solution to the 50-mL line marked on the Wash 3 Bottle, then cap the bottle.

## Begin the initialization

**Note:**

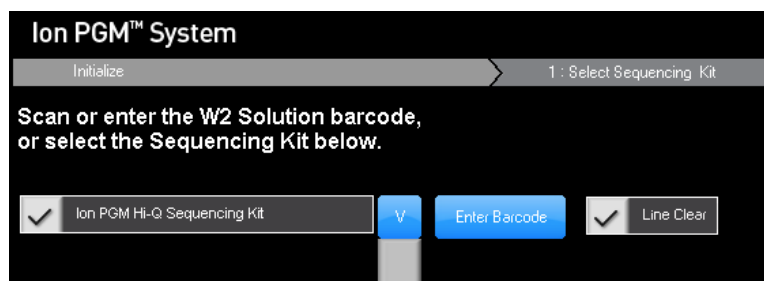
- Do not remove the old sipper tubes from the dNTP ports until instructed to do so.
- Load the bottles as quickly as possible to prevent atmospheric CO<sub>2</sub> from reducing the pH of the Wash 2 Bottle solution.

---

**IMPORTANT!** Do not let the new sipper tubes touch any surfaces.

---

- On the main menu, press **Initialize**.
- Make the following selections in the next screen, then press **Next**:
  - Click **Enter barcode** to scan or enter the barcode on the Ion PGM™ Hi-Q™ Sequencing W2 Solution bottle, or the 2D barcode on the Ion PGM™ Sequencing Solutions box.
  - Alternatively, select the checkbox for the **Ion PGM™ Hi-Q™ Sequencing Kit** from the dropdown list.
  - In the same screen, if you routinely experience clogging during initialization, select the **Line Clear** checkbox to clear any blockage in the fluid lines before initialization. This is optional.



**IMPORTANT!** Be careful to select the correct kit, to ensure proper pH adjustment.

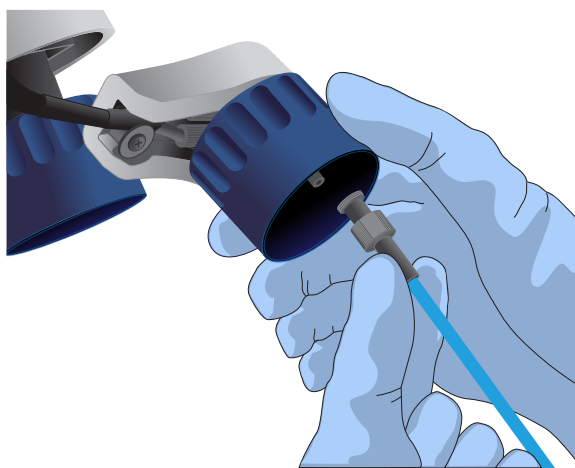
After you press **Next**, the system will check the gas pressure.

3. Following the gas pressure check:

Result	Action
<b>If the pressure is sufficient</b>	Ensure that the cleaning chip, reagent sipper tubes, and collection trays are in place, and press <b>Next</b> to start the initialization.
<b>If the pressure is low</b>	Press <b>Yes</b> to re-check the pressure. If the pressure remains low, contact Technical Support.

4. Wearing clean gloves, firmly attach a new, long gray sipper to the cap in the W2 position.

**IMPORTANT!** Do not let the sipper touch any surfaces, and firmly attach the sipper to the port. Loosely attached sippers can adversely affect results.



5. Immediately attach the prepared Wash 2 Bottle in the W2 position, then tighten the cap. Press **Next**.
6. Change gloves and firmly install new sipper tubes (short gray) in the caps in the W1 and W3 positions.
7. Immediately attach the prepared Wash 1 and 3 Bottles, then tighten the caps. Press **Next**.

8. If you selected the **Line Clear** checkbox in the earlier screen, press **Next**, then follow the touchscreen prompts to perform the line clear procedure. At the beginning and end of the procedure, you are prompted to select one of the following:

Option	Description
<b>Press Line Clear</b>	To start a new line clear procedure
<b>Press Re-flow</b>	To retest the lines after you have performed a line clear
<b>Press Auto pH</b>	If the lines are clear and you are ready to continue with initialization

9. Following line clear, or if you did not select that option, the sequencer begins adjusting the pH of the W2 Solution, which takes ~30 minutes. After 15 minutes, check the instrument touchscreen to confirm that initialization is proceeding normally.

**Note:**

- If an error occurs during the automatic pH process, note the error message and proceed to “Initialization—Auto pH errors” on page 94.
- During the process, you can start preparing the Reagent Bottles with dNTPs as described in the next section.

## Prepare the 50-mL Reagent Bottles with dNTPs

1. Use the labels provided with the kit to label four new Reagent Bottles as dGTP, dCTP, dATP, and dTTP.
2. Confirm that no ice crystals are visible in each thawed dNTP stock solution. Vortex each tube to mix, and centrifuge to collect the contents. Keep the dNTP stock solutions on ice throughout this procedure.

---

**IMPORTANT!** To avoid cross-contamination in the next step, open only one dNTP stock tube at a time and use a fresh pipette tip for each aliquot.

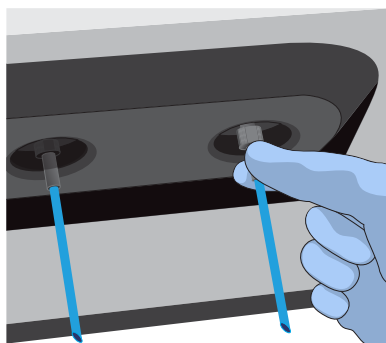
---

3. Using separate filtered pipette tips and clean gloves, carefully transfer 20 µL of each dNTP stock solution into its respective Reagent Bottle.
4. Cap each Reagent Bottle and store on ice until you are ready to attach it to the instrument. Place the remaining dNTP stocks back into –20°C for storage.

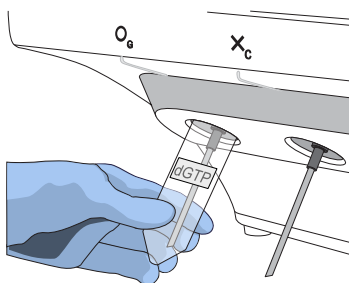
## Attach the sipper tubes and Reagent Bottles

1. After the wash solutions have initialized, follow the touchscreen prompts to remove the used sipper tubes and collection trays from the dNTP ports.
2. Change gloves, then firmly insert a new sipper tube (blue) into each dNTP port. Do not let the sipper touch any surfaces.

**IMPORTANT!** Be careful to firmly push each sipper onto the port. Loosely attached sippers may adversely affect results.



3. Attach each prepared Reagent Bottle to the correct dNTP port (e.g., the dGTP tube on the port marked "G") and tighten firmly by hand until snug. Press **Next**.



**Note:** The instrument checks the pressure of the Reagent Bottles and Wash Bottles. If a bottle leaks, check that it is tightly attached to the instrument. If it continues to leak, replace it. If the instrument still does not pass the leak check, contact Technical Support.

4. Follow the touchscreen prompts to complete initialization. The instrument will fill each Reagent Bottle with 40 mL of W2 Solution.
5. At the end of initialization, Ion PGM™ System will measure the pH of the reagents:
  - If every reagent is in the target pH range, a green **Passed** screen will be displayed.
  - If a red failure screen appears, see Appendix A, "Troubleshooting".
6. Press **Next** to finish the initialization process and return to the main menu.
7. Proceed to the appropriate sequencing protocol for your chip type.



# Load the chip and start the sequencing run

■ Materials required .....	64
■ Guidelines for sequencing .....	65
■ Before you begin .....	65
■ Optional: Prepare Ion Sphere™ Test Fragments .....	65
■ Add controls to the enriched, template-positive ISPs .....	65
■ Anneal the Sequencing Primer .....	66
■ Perform Chip Check .....	67
■ Bind the Sequencing Polymerase to the ISPs .....	69
■ Prepare and load the chip .....	70
■ Select the Planned Run and perform the run .....	74

Use the following chip loading and sequencing protocol for all Ion PGM™ chip types.

## Materials required

### Materials provided in the kit

- Sequencing Primer
- Control Ion Sphere™ Particles
- Annealing Buffer
- Ion PGM™ Hi-Q™ Sequencing Polymerase

### Other materials and equipment

- Ion Chip kit: Ion 318™ Chip v2 BC, Ion 316™ Chip v2 BC, or Ion 314™ Chip v2 BC
- Enriched template-positive ISPs
- 0.2-mL PCR tube (non-polystyrene)
- Rainin™ SR-L200F pipette and tips
- Vortex mixer
- Ion Chip™ Minifuge
- Thermal cycler with heated lid (programmed at 95°C for 2 minutes and 37°C for 2 minutes)
- Barcode scanner (included with the Ion PGM™ System)



## Guidelines for sequencing

- Ion PGM™ Sequencer: When performing multiple sequencing runs in the same instrument initialization, the first run should be started within 1 hour after initialization, and the last run must be started within 24 hours after initialization.
- Ion PGM™ Sequencer: If you press the **Abort** button on the touchscreen during any procedure, the touchscreen can freeze, requiring that you restart the sequencer.
- The ISPs are difficult to see. To avoid aspirating the particles:
  - When centrifuging the ISPs, orient the tab of the tube lid so that it is pointing away from the center of the centrifuge, to indicate where the pellet will be formed.
  - Always remove supernatant from the tube from the top down.

## Before you begin

1. Thaw the Sequencing Primer on ice.
2. Update the Torrent Suite™ System and Ion PGM™ System software to the latest version available, if needed.

## Optional: Prepare Ion Sphere™ Test Fragments

If you are performing an installation or troubleshooting sequencing run:

1. Vortex the Ion Sphere™ Test Fragments from the Ion PGM™ Controls Kit (Cat. No. 4480449), then pulse-centrifuge in a microcentrifuge for 2 seconds before taking aliquots.
2. Add 5 µL of Ion Sphere™ Test Fragments to 100 µL of Annealing Buffer in a 0.2-mL non-polystyrene PCR tube.

Skip directly to "Anneal the Sequencing Primer".

## Add controls to the enriched, template-positive ISPs

1. Vortex the Control Ion Sphere™ Particles, then pulse-centrifuge in a picofuge for 2 seconds before taking aliquots.
2. Add 5 µL of Control ISPs directly to the entire volume of enriched, template-positive ISPs (prepared using your template preparation method) in a 0.2-mL non-polystyrene PCR tube.

Proceed to "Anneal the Sequencing Primer".

## Anneal the Sequencing Primer

1. Mix the tube containing the ISPs (or test fragments) by thoroughly pipetting up and down.
  2. Place the tube in a microcentrifuge with an appropriate tube adapter. Orient the tab of the tube lid so that it is pointing away from the center of the centrifuge, to indicate where the pellet will be formed.
  3. Centrifuge for 3 minutes at  $14,000 \times g$ .
  4. Keeping the pipette plunger depressed, insert a pipette tip into the tube containing the pelleted ISPs and carefully remove the supernatant from the top down, avoiding the side of the tube with the pellet (i.e., the side with the tab on the tube lid). Discard the supernatant. Leave  $\sim 15 \mu\text{L}$  in the tube (visually compare to  $15 \mu\text{L}$  of liquid in a separate tube).
  5. Ensure that the Sequencing Primer is completely thawed prior to use (no ice crystals should be visible).
  6. Vortex the primer for 5 seconds, then pulse spin in a microfuge for 3–5 seconds to collect the contents. Leave on ice until ready to use.
  7. Add  $12 \mu\text{L}$  of Sequencing Primer to the ISPs, and confirm that the total volume is  $27 \mu\text{L}$  (add Annealing Buffer if necessary).
  8. Pipet the mixture up and down thoroughly to disrupt the pellet.
- 
- IMPORTANT!** Make sure that the pipette tip is at the bottom of the tube during mixing to avoid introducing air bubbles into the sample.
- 
9. Program a thermal cycler for  $95^{\circ}\text{C}$  for 2 minutes and then  $37^{\circ}\text{C}$  for 2 minutes, using the heated lid option.
  10. Place the tube in the thermal cycler and run the program. After cycling, the reaction can remain in the cycler at room temperature ( $20\text{--}30^{\circ}\text{C}$ ) while you proceed with Chip Check.

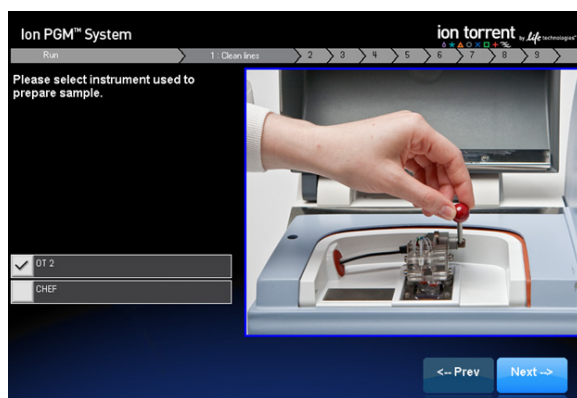
## Perform Chip Check

Chip Check tests the chip and ensures that it is functioning properly before loading the sample.

### IMPORTANT!

- To avoid damage due to electrostatic discharge (ESD), **do not place the chip directly on the bench or any other surface**. Always place the chip either on the grounding plate on the Ion PGM™ Sequencer or in the Ion Chip™ Minifuge adapter bucket.
- To avoid ESD damage, **do not wear gloves** when transferring chips on and off the instrument.

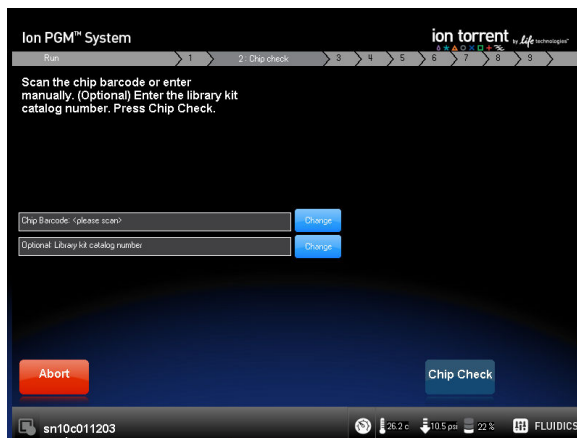
1. On the main menu of the Ion PGM™ Sequencer touchscreen, press **Run**. Remove the waste bottle and completely empty it. Press **Next**.
2. When prompted to insert a cleaning chip, use the same used chip that was used for initialization. Press **Next** to clean the fluid lines.
3. When prompted, select the instrument that you used to prepare the template-positive ISPs. Then press **Next**.



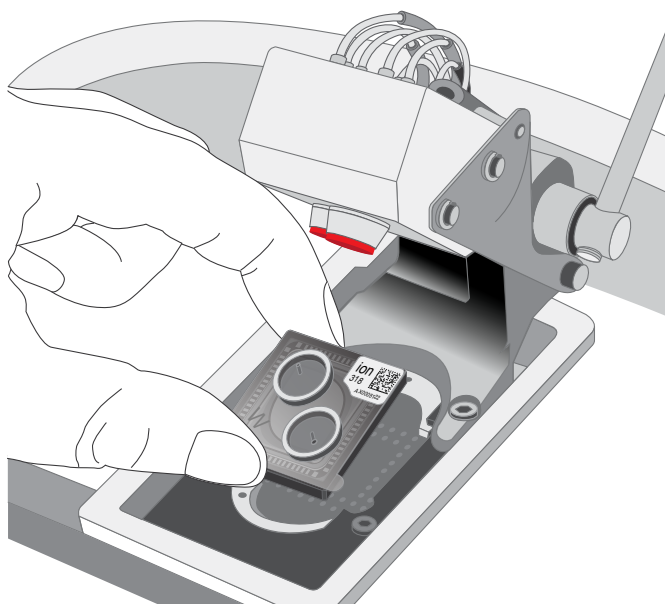
4. Remove gloves, then ground yourself by touching the grounding pad on the sequencer. Remove a new chip from its packaging, then label it to identify the experiment (save the chip package). Press **Next**.

- When prompted, use the scanner to scan the barcode located on the new chip, or press **Change** to enter the barcode manually. Optionally, you can also enter the library kit catalog number.

**Note:** A chip cannot be run without scanning or entering the barcode.



- Replace the old chip in the chip socket with the new one. Close the chip clamp, then press **Next**.



- Press **Chip Check**. During the initial part of Chip Check, visually inspect the chip in the clamp for leaks.

**Note:**

- If there is a leak, press the **Abort** button immediately to stop the flow to the chip. Proceed to Appendix A, "Troubleshooting".
- The chip socket can be damaged by rubbing or wiping its surface.** Never rub or wipe the socket to clean up leaks. See Appendix A, "Troubleshooting" for more information.

8. When Chip Check is complete:

- If the chip passes, press **Next**.
- If the chip fails, open the chip clamp, reseal the chip in the socket, close the clamp, and press **Calibrate** to repeat the procedure. If the chip passes, press **Next**. If the chip still fails, press **Main Menu** and restart the experiment with a new chip. See Appendix A, "Troubleshooting" for more information.

**Note:** To return *damaged* chips, contact Technical Support.

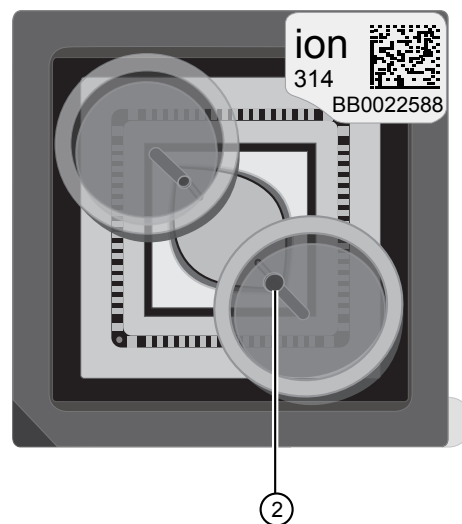
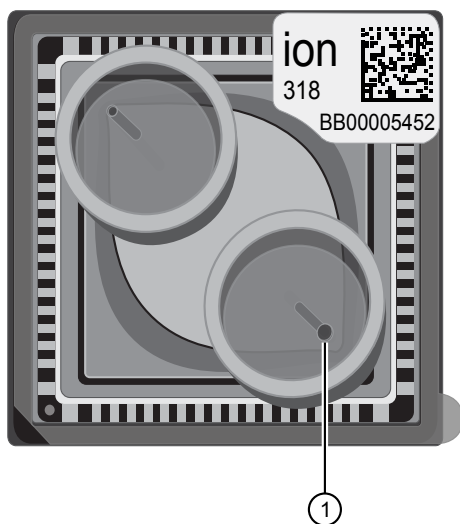
9. Following a successful Chip Check, empty the waste bottle, then select the **Waste bottle is empty** checkbox on the touchscreen. Press **Next**.

## Bind the Sequencing Polymerase to the ISPs

1. Remove the Ion PGM™ Hi-Q™ Sequencing Polymerase from storage and flick mix with your finger tip 4 times. Pulse-centrifuge for 3–5 seconds. Place on ice.
2. After annealing the Sequencing Primer, remove the ISPs from the thermal cycler, then add 3 µL of Ion PGM™ Hi-Q™ Sequencing Polymerase to the ISPs, for a total final volume of 30 µL.
3. Pipet the sample up and down to mix, then incubate at room temperature for 5 minutes.

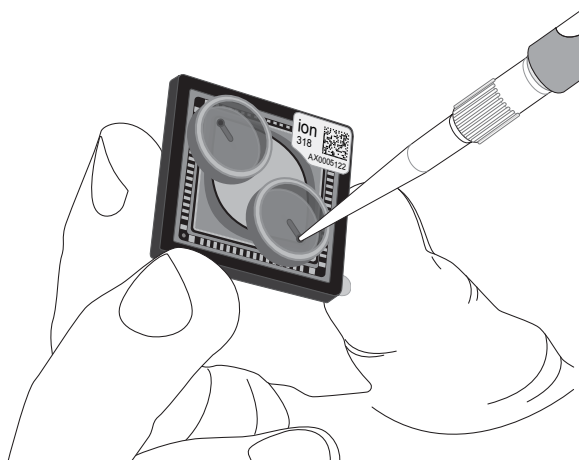
## Prepare and load the chip

Remove liquid  
from the chip



- ① Ion 318™ Chip v2 BC/Ion 316™ Chip v2 BC      ② Ion 314™ Chip v2 BC loading port

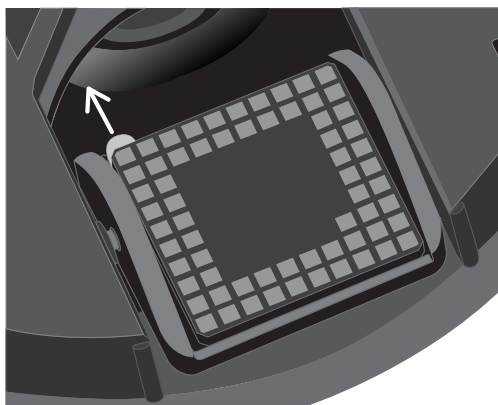
1. Following chip calibration, remove the new chip from the Ion PGM™ Sequencer. Insert a used chip in the chip clamp while loading the new chip.
2. Tilt the new chip at a 45° angle so that the loading port is the lower port.



3. Insert the pipette tip firmly into the loading port, then remove as much liquid as possible from the loading port. Discard the liquid.

**IMPORTANT!** For the next steps, if you are preparing one chip at a time, balance the chip in the Ion Chip™ Minifuge rotor with a used chip of the same chip type and orientation. Be careful to balance an upside-down chip with another upside-down chip. Mark the used chip with a laboratory marker to differentiate it from the new chip.

4. Place the chip **upside-down** in the minifuge bucket, then transfer the bucket **with the chip tab pointing in** (toward the center of the minifuge). Balance the bucket with another chip.



5. Centrifuge for 5 seconds to empty the chip completely.



**CAUTION!** Allow the minifuge to come to a complete stop before opening the lid.

6. Remove the chip from the bucket, then wipe the bucket with a disposable wipe to remove any liquid. Place the chip right-side up in the bucket.

## Load the chip

1. Place the chip in the bucket on a firm, flat surface. Following polymerase incubation, load the chip with following volume of prepared ISPs using the listed pipettes, or equivalent, depending on your chip type. We recommend using a P20 pipette for Ion 314™ Chips for optimal loading.

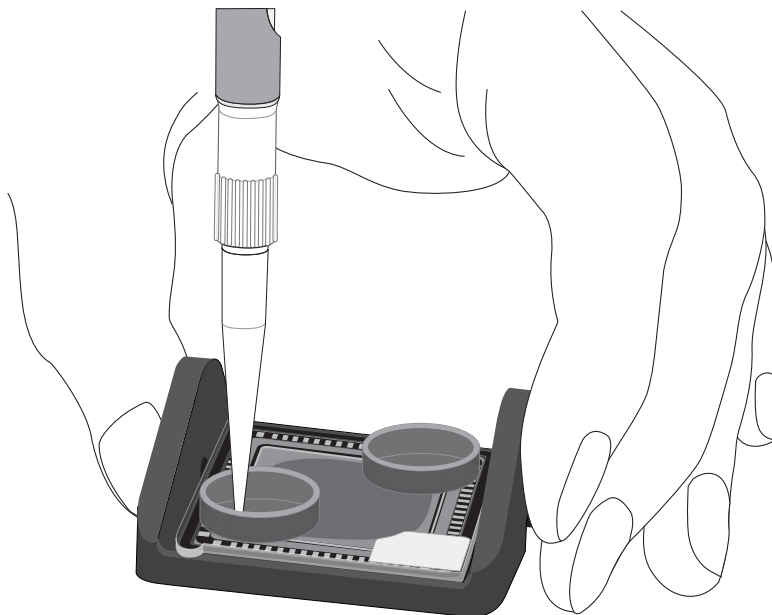
Chip	Volume to load	Recommended pipette <sup>[1]</sup>
Ion 316™ or Ion 318™ Chip	Entire volume (~30 µL)	Rainin™ Pipet-Lite™ LTS L-100XLS, 10–100 µL
Ion 314™ Chip	10 µL	Rainin™ Pipet-Lite™ LTS L-20XLS, 2–20 µL

<sup>[1]</sup> Alternatives from Gilson and Eppendorf can be used.

**Note:** For Ion 314™ Chips, the remaining volume of ISPs can be used to prepare another chip, provided the chip can be loaded and sequenced in parallel with the first chip. After polymerase incubation, ISPs should be loaded and sequencing initiated within ~15 minutes.

2. Insert the tip firmly into the loading port of the chip.

**Note:** When loading the ISPs into the chip, keep the pipette tip at a 90° angle to the chip, press the tip firmly into the circular loading port, and apply gentle pressure between the pipette tip and chip.



3. With the pipette unlocked, apply gentle pressure between the tip and chip and slowly dial down the pipette (~1  $\mu\text{L}$  per second) to deposit the ISPs. To avoid introducing bubbles into the chip, leave a small amount in the pipette tip (~0.5  $\mu\text{L}$ ).

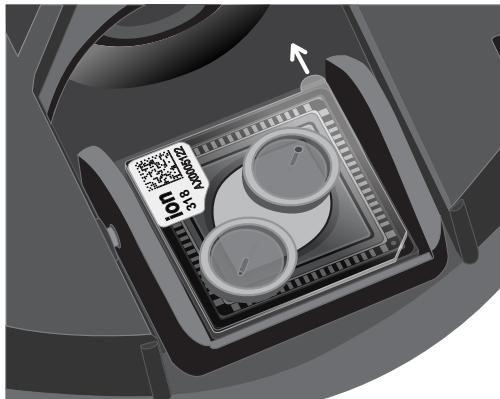


**Note:** Do not remove the pipette tip from the port during the dial-down process, because removal can introduce air bubbles and inhibit loading.

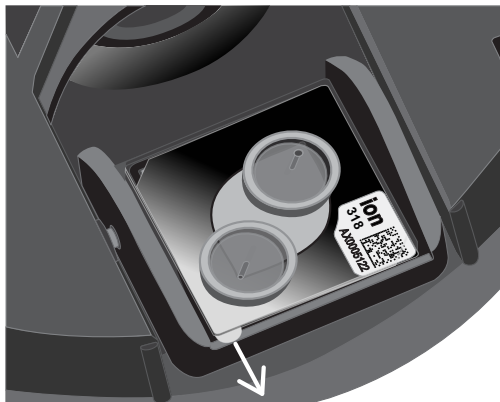
4. Remove, then discard any displaced liquid from the other port of the chip.



5. Transfer the chip in the bucket to the minifuge with the chip tab **pointing in** (toward the center of the minifuge), then centrifuge for 30 seconds.



6. Turn the chip so that the chip tab is **pointing out** (away from the center of the minifuge), then centrifuge for 30 seconds.



7. Remove the bucket from the minifuge, then place it on a flat surface. Set the volume of the pipettor as follows, depending on your chip type:
  - **Ion 316™ or Ion 318™ Chip:** 25  $\mu$ L
  - **Ion 314™ Chip:** 5  $\mu$ L
8. Tilt the chip at a 45° angle so that the loading port is the lower port, then insert the pipette tip into the loading port.
9. Without removing the tip, slowly pipet the sample out and then back into the chip one time. **Pipet slowly to avoid creating bubbles.**
10. Slowly remove as much liquid as possible from the chip by dialing the pipette. Discard the liquid.
11. Turn the chip upside-down in the bucket, transfer it back to the minifuge, then centrifuge upside-down for 5 seconds. Remove and discard any liquid.

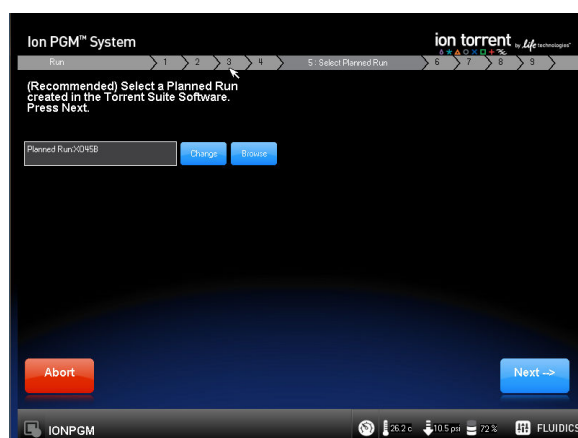
12. If some liquid remains in the chip, lightly and rapidly tap the point of the chip tab against the benchtop a few times, then remove and discard any collected liquid. Do not flush the chip.
13. When chip loading is complete, press **Next** on the touchscreen, then proceed immediately to performing the run.

## Select the Planned Run and perform the run

### Select the Planned Run

1. Press **Browse** next to the **Planned Run** field and select the name of the plan you created, then touch **Next**.

**Note:** The Ion PGM™ Sequencer automatically populates this field for barcoded Ion chips.



2. Confirm that the settings are correct. If necessary, make any changes using the touchscreen controls.



**Note:** If the number of flows (cycles) to be run cannot be selected, there may not be enough disk space to store the experiment data. Touch **Data Mngt** to start the Data Management application (this can also be accessed from the Tools Menu) and delete old runs from the Ion PGM™ System.

## Perform the run

1. After you enter the Planned Run, press **Next** to verify the experimental setup. Press **OK** to confirm the settings or press **Cancel** to return to the touchscreen to adjust the settings.
2. When prompted by the instrument, load and clamp the chip, then press **Next**.
3. At the beginning of the run, visually inspect the chip in the clamp for leaks before closing the cover. The instrument will flush any loose ISPs from the chip and begin calibrating the chip.
4. When the calibration is complete (~1 minute), the touchscreen will indicate whether calibration was successful.
  - If the chip passes calibration, press **Next** to proceed with the sequencing run.
  - If the chip fails calibration, see “Error message: Calibration FAILED” on page 93.
5. After 60 seconds, the run will automatically begin, or press **Next** to begin the run immediately.

---

**IMPORTANT!** During a run, avoid touching the instrument and any of the attached bottles or tubes, as this may reduce the quality of the measurements.

---

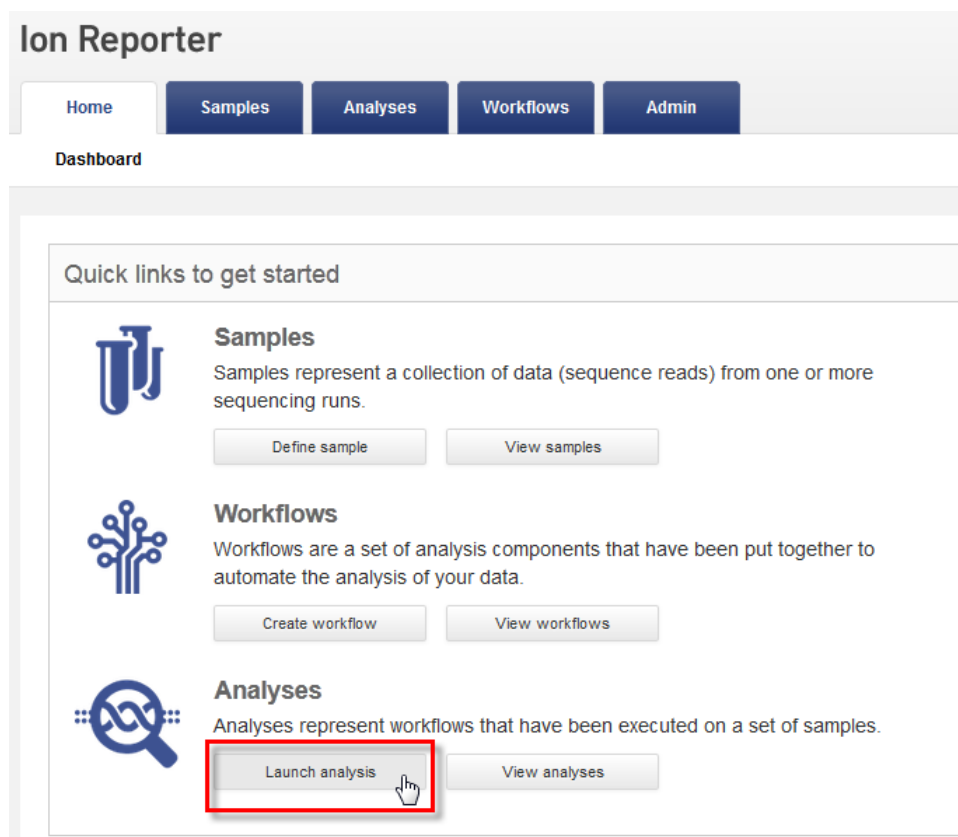
6. When the run is complete, leave the chip in place, then touch **Next** to return to the Main Menu. You can then remove the chip and proceed with another run or perform a cleaning/initializing if required.

**Note:** See “Cleaning schedule” on page 55 to determine whether cleaning is required after the run.

## Launch an Ion Reporter™ analysis

If you are analyzing your samples manually, follow the instructions below. If you planned your run for automatic analysis with Ion Reporter™ Software, proceed to step 6.

1. Import your samples into Ion Reporter™ Software using the Ion Reporter™ Uploader plugin.  
See the *Ion Reporter™ Software Help Guide* (Pub. No. MAN0013516) for detailed instructions for importing your samples and defining them in Ion Reporter™ Software.
2. In the Ion Reporter™ **Home** tab, click **Launch analysis** after the Ion Reporter™ Uploader plugin has completed.



3. In the **Workflow** tab, select **Aneuploidy** from the **Applications** menu, and **Ion** from the **Workflow** menu to view the Ion aneuploidy workflows. Select **ReproSeq Low-pass whole-genome aneuploidy** from the list, then click **Next**.

#### Launch Analysis

Workflow Samples Plugins

Select the workflow you wish to launch. [Learn more...](#)

Aneuploidy Workflow Target Group 5.2 Reference Search

ion	Application	Workflow Name	Version	Reference	Sample Group	Modified On
ion	Aneuploidy	ReproSeq Low-pass whole-genome aneuploidy	5.2	hg19	Single	Sep 27 2016 03:13 AM
ion	Aneuploidy	Low-pass whole-genome aneuploidy	5.2	hg19	Single	Sep 27 2016 03:13 AM

1 - 2 of 2 items

Cancel Next

**Note:** The ReproSeq Low-Coverage Whole-Genome Baseline is auto-loaded in the ReproSeq Low-pass whole-genome aneuploidy workflow, making creation of a custom baseline and workflow for your Ion ReproSeq™ analyses unnecessary.

4. In the **Samples** tab, select the sample(s) you want to run in your analysis from the list.

#### Launch Analysis

Workflow Samples Plugins

Select the sample you wish to analyze. You can select multiple samples and each one will be treated as a separate analysis. [Learn more...](#)

Samples Search

	Sample	Gender	Sample Type	Role	Imported By	Imported On
<input checked="" type="checkbox"/>	WSSample_1493a0d4-1d2f-466a-a402-9b5835d0046a	Male	Unknown	proband	User, Ion	Oct 24 2016 05:04 AM
<input checked="" type="checkbox"/>	WSSample_66521b70-dae9-4143-849b-7b12d4344586	Male	Unknown	proband	User, Ion	Oct 24 2016 05:03 AM
<input type="checkbox"/>	cv_bam_IRUCLL_DNA_1	Unknown	Unknown	proband	User, Ion	Oct 24 2016 01:41 AM

5. Click **Next** twice to advance through the **Plugins** tab to **Confirm & Launch**.

6. In the **Confirm & Launch** tab, enter a name for your analysis if you want to change the default name, then click **Launch Analysis**.

**Launch Analysis**

Workflow Samples Plugins **Confirm & Launch**

**Analysis ready to launch!**

Review the selected options, name your analysis and then launch it.

Analysis Name: WSSample\_1493a0d4-1d2f-466a-a402-9b5835d0046a\_c949\_2016-10-24-16-52-9t  
(WSSample\_1493a0d4-1d2f-466a-a402-9b5835d0046a)

Description: Optional

**Launch Analysis**

← Previous Cancel

**Summary**

Application: Aneuploidy

Workflow: ReproSeq Low-pass whole-genome aneuploidy

Annotations: Aneuploidy

Filters: CNVs of Confidence >= 0.1 - Germline - CNVs only

Samples: 1 Sample

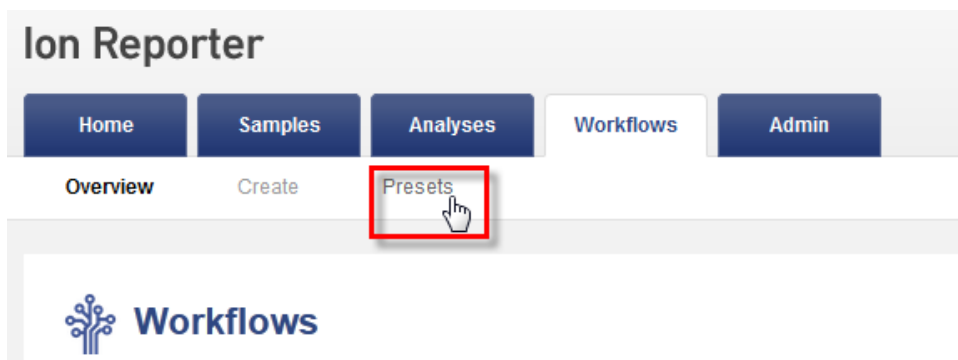
Price: \$0.00 USD

7. Review your results by selecting from the Analyses list after navigating to **Analysis ▶ Overview**. See the *Ion Reporter™ Software Help Guide* (Pub. No. MAN0013516) for further details.

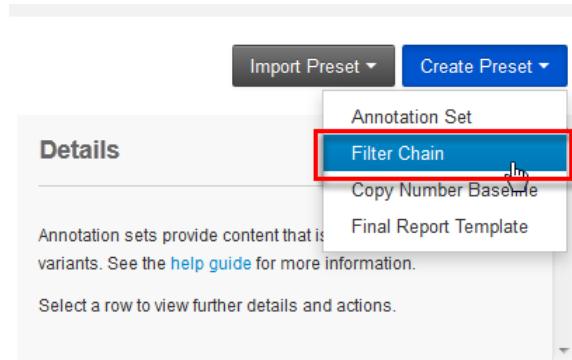
## Edit the Ion Reporter™ ReproSeq Low-pass whole-genome aneuploidy workflow

The ReproSeq Low-pass whole-genome aneuploidy workflow has the CNV Confidence Range set at 0.1 to 1.0E7 as default. To specify another range for more stringent (1.0 to 1.0E7) or less stringent (0.01 to 1.0E7) CNV filtering, first create a new Filter Chain, then make a copy of the workflow and edit it to add the new Filter Chain.

1. In the Ion Reporter™ **Workflows** tab, click **Presets**.



- In the **Create Preset** dropdown list, select **Filter Chain**.



- In the dialog, select **CNV Confidence Range - CNVs Only** from the dropdown list on the left, enter the Confidence Range, then click **Set**.

### Create Filter Chain

Name: Required

Description: Optional

Reference: GRCh38, hg19

CNV Confidence Range - CNVs

Range: 0.0 ← → 1.0E7

From: 1.0

To: 1.0E7

☒ Include boundary values

☐ Include unannotated variants

**Set**

- Enter a name for the Filter Chain, then click **Save**.

### Create Filter Chain

Name: ReproSeq\_10142016

Description: Optional

Reference: GRCh38, hg19

CNV Confidence Range - CNVs

Range: 0.0 ← → 1.0E7

From: 1.0

To: 1.0E7

☒ Include boundary values

☐ Include unannotated variants

**Set**

FilterChain Query: CNV Confidence Range - CNVs Only

Selected Filters:

Name	Value
CNV Confidence Range - CNVs Only	1.0 ≤ CNV Confidence Range - CNVs Only ≤ 1.0E7

**Save**

5. Select the **ReproSeq Low-pass whole-genome aneuploidy** workflow from the list of workflows in the Ion Reporter™ **Workflows** tab, then click **Copy** in the **Actions**▼ dropdown list.

Application	Workflow Name	Version	Reference	Sample Group	Modified On
Ion	ReproSeq Low-pass whole-genome aneuploidy	5.2	hg19	Single	Oct 23 2016 10:42 PM
Ion	Low-pass whole-genome aneuploidy	5.2	hg19	Single	Oct 23 2016 10:42 PM

Selected Workflows: 2

Details: Copy, Launch Analysis, Tag for IRU

6. In the Filters tab of the Edit Workflow wizard, select your Filter Chain from the dropdown list, then click **Next**.

**Edit Workflow** copy of ReproSeq Low-pass whole-genome aneuploidy\_07c6l

Application Reference Annotation **Filters** Copy Number Plugins Final Report

Filters allow you to automatically remove variants that meet certain criteria. Select the filter set you wish to apply to the detected variants. [Learn more...](#)

**Filter Chains**

- Ion CNVs of Confidence >= 0.0...
- Ion OncoPrint Variants
- Ion OncoPrint Variants, 5% CI CNV ploidy => gain of 2 over normal
- phyloP\_filter\_test
- ReproSeq\_10142016**

Next

7. Advance through the remaining tabs of the wizard to make any additional edits, then name your edited workflow in the **Confirm** tab, then click **Save Workflow**.

**Edit Workflow** copy of ReproSeq Low-pass whole-genome aneuploidy\_07c60056-6cba-40f3-9d3f-ec5472888951

Application Reference Annotation Filters Copy Number Plugins Final Report Parameters **Confirm**

**Workflow configured!**

Review the selected options, name your workflow and then save it.

Workflow Name: ReproSeq Low-pass whole-genome aneuploidy\_edit1

Description: Detects aneuploidies and large chromosome abnormalities from a single whole-genome sample with low coverage (minimum 0.01x). Normalization is done using...

**Save Workflow**

Previous Cancel

**Summary**

- Application: Aneuploidy
- Sample Group: Single
- Reference: hg19
- Annotations: Aneuploidy
- Filters: ReproSeq\_10142016
- Copy Number: ReproSeq Low-Coverage Whole-Genome Baseline
- Report Template: Default Final Report Template

**Note:** See the *Ion Reporter™ Software Help Guide* (Pub. No. MAN0013516) for further details.



## Visualize results with IGV Light viewer

You can visualize aneuploidy results and generate reports with a streamlined IGV Light viewer. See the *Ion Reporter™ Software Help Guide* for further details.

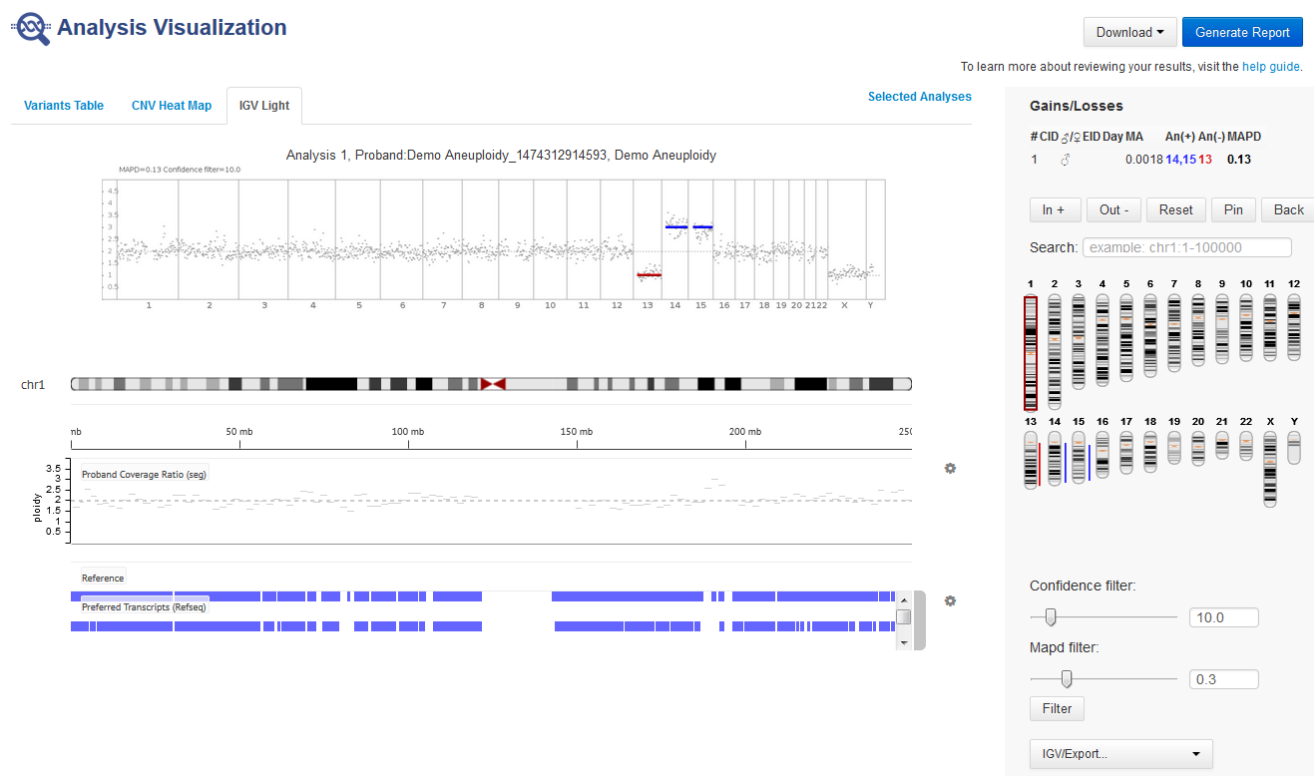
1. On the Ion Reporter™ home page, navigate to **Analyses ▶ Overview**.
2. Select one or more aneuploidy analyses, then click **Visualize** from the **Actions** dropdown list.

The screenshot shows the 'Analyses' overview page in Ion Reporter. At the top, there are filters for Status, Aneuploidy, Version (5.2), and Reference. Below these is a table of analyses. The first row is selected, and its 'Actions' dropdown menu is open, showing options like Archive, Audit Log, Delete, Download Filtered Variants, Download All Variants, Edit Analysis Results, Download Logs, Send to Report Role, Share, View QC Report, and Visualize. The 'Visualize' option is highlighted with a red box.

	Analysis	Version	Reference	Stage	Project	Workflow	Created On	Status
<input checked="" type="checkbox"/>	IRUCLI_Automation_05-10-2016-11-50-38-550_Aneuploidy_3c4f459b-2a4b-4822-b82e-83ca0de19381	5.2	hg19	Interpretation Assignment	Unknown	ReproSeq Low-pass whole-genome aneuploidy	Oct 05 2016 11:54 AM	Successful
<input type="checkbox"/>	Automation_BAC_05-10-2016-11-40-52-22	5.2	hg19	Interpretation Assignment		ReproSeq Low-pass whole-genome aneuploidy	Oct 05 2016 11:41 AM	Successful
<input type="checkbox"/>	SampleDefineKVJe_c1149_2016-10-05-04-18-483	5.2	hg19	Interpretation Assignment		PS_Cus_ReproSeq_aneuploidy_rmDup_T rue_12Sep2016	Oct 05 2016 04:19 AM	Successful
<input type="checkbox"/>	Automation_SampleDefine_05-10-2016-04-10-02-826_c1155_2016-10-05-04-10-644	5.2	hg19	Analysis		PS_Cus_ReproSeq_aneuploidy_rmDup_T rue_12Sep2016	Oct 05 2016 04:11 AM	Aborted
<input type="checkbox"/>	Demo Aneuploidy_1474312916203	5.2	hg19	Interpretation Assignment		ReproSeq Low-pass whole-genome aneuploidy	Sep 19 2016 12:22 PM	Successful

**Note:** You can also click **Visualize** (to the left of the **Launch Analysis** button) after selecting two or more analyses.

3. Click the **IGV Light** tab. A copy number histogram for each analysis selected appears, along with ploidy maps for selected chromosomes or chromosome regions, and karyograms showing copy number gains and losses.



**Note:** If you select multiple analyses generated from workflows with different confidence filter settings, the software will use the lowest confidence filter setting among the analyses and apply it across the group.

4. Select **Open in IGV (download .jnlp file)** from the **IGV/Export...** dropdown list, or click **Download** at the upper right to download a file from which to launch the IGV application. You can also click **Generate Report** to produce a full report with graphics that can be downloaded as a pdf.

To learn more about reviewing your results, visit the [help guide](#).

**Gains/Losses**

#	CID	EID	Day	MA	An(+)	An(-)	MAPD
1					0.0018	14,15,13	0.135

In + Out - Reset Pin Back

Search:

1 2 3 4 5 6 7 8 9 10 11 12  
13 14 15 16 17 18 19 20 21 22 X Y

Confidence filter:

Mapd filter:

Filter

IGV/Export...  
 Open in IGV (download .jnlp file)  
 Export whole genome view (if available) only to PNG  
 Export IGV tracks only to PNG  
 Show IGV preferences



# Troubleshooting

## Ion SingleSeq™ library preparation

Observation	Possible cause	Recommended action
Library yield is low.	Inefficient amplification of genomic DNA.	Use the Human CEPH Genomic DNA Control at 15–60 pg input as a positive control in pre-amplification and amplification reactions to determine whether DNA is amplifiable in your system.
		Use real-time PCR to monitor amplification.
	Insufficient starting material due to loss or mishandling of cells.	Increase the number of cells in your sample.
		Keep the cells on ice.
	Over-drying of AMPure™ XP Reagent during cleanup.	Do not insert pipette tip into sample – cell(s) may adhere to the tip after addition of reagent.
Barcode balance in a pool is poor.	Variation in gDNA input in a pool of samples is too high.	Do not dry the AMPure™ XP beads more than 4 minutes.
		Avoid pooling single cell samples with multiple cell samples; i.e. pool single-cell samples together and pool multiple-cell samples in a second pool if possible.




## Ion OneTouch™ ES

For Ion OneTouch™ ES vertical and horizontal axis calibration and residual volume test procedures, see Chapter 3 of the *Ion OneTouch™ 2 System User Guide* (Pub. No. MAN0014388).

Observation	Possible cause	Recommended action
Excessive foaming occurs	<ul style="list-style-type: none"><li>Instrument is improperly calibrated resulting in inadequate volume in one or more wells.</li><li>Fitting is loose.</li><li>Pipette tip is cracked.</li></ul>	<ol style="list-style-type: none"><li>Use the recommended volumes for all wells.</li><li>Ensure that fittings are tight, especially at the elbow fitting, and the pipette tip is not cracked.</li><li>If necessary, perform the residual volume test. If the residual volume test fails, then calibrate the instrument.</li></ol>



Observation	Possible cause	Recommended action
Brown pellet is present in centrifuged tube of enriched ISPs	Residual Dynabeads™ MyOne™ Streptavidin C1 Beads are present.	<ol style="list-style-type: none"> <li>1. Pipet the suspension with the brown pellet up and down 10 times to resuspend the pellet.</li> <li>2. Place the 0.2-mL PCR tube against a magnet such as a DynaMag™-2 magnet for 4 minutes.</li> <li>3. Transfer the supernatant with the enriched ISPs to a new 0.2-mL PCR tube without disturbing the pellet of Dynabeads™ MyOne™ Streptavidin C1 Beads.</li> <li>4. Sequence or store the enriched ISPs.</li> </ol>

Observation	Possible cause	Recommended action
E12, E22, or E23 errors display during the run or during calibration	Calibration values are out of range.	<ol style="list-style-type: none"> <li>1. Power OFF the instrument and wait 3 seconds.</li> <li>2. While holding down <b>Vert. Adjust</b>, power ON the instrument. This step restores the factory default settings.</li> <li>3. Recalibrate the vertical axis: <p><b>Note:</b> The default setting for the vertical axis is 310. If the setting is &lt;310, the instrument will likely display an error, because the Tip Arm position is too high.</p> <ol style="list-style-type: none"> <li>a. Press the  (minus) button to lower the Tip Arm until the tip touches the shelf.</li> <li>b. Press the  (minus) button 8 more times. Typical vertical axis settings are ~340–370.</li> </ol> </li> <li>4. Recalibrate the horizontal axis: Press the  (plus) button to move the Tip Arm to the right until the tip touches the left tab of the strip. <p><b>Note:</b> The default setting for the horizontal axis is 625. Typical horizontal axis settings are ~640–670.</p> </li> </ol>
	AC line voltage module is installed incorrectly.	<ol style="list-style-type: none"> <li>1. Determine the voltage of the electrical outlet to plug in the Ion OneTouch™ ES.</li> <li>2. Align the arrow by the correct voltage on the AC line voltage module with the adjacent white arrow in the lower-right corner of the fuse socket.</li> </ol> <p>If the AC line voltage module is installed incorrectly:</p> <ol style="list-style-type: none"> <li>1. Gently remove the module with your fingernail or a small flathead screwdriver.</li> <li>2. Rotate the module so that the correct voltage on the module is aligned and adjacent to the white arrow in the lower right-hand corner of the fuse socket.</li> <li>3. Insert the AC line voltage module into the fuse socket.</li> </ol>



Observation	Possible cause	Recommended action
E12 or E22 error is displayed when the unit is initializing	<ul style="list-style-type: none"> <li>Fuse is installed incorrectly.</li> <li>Unit is below operating temperature.</li> <li>Program or calibration setting is bad, <i>or</i></li> <li>Tip Arm is not moving.</li> </ul>	<ol style="list-style-type: none"> <li>Ensure that the fuse module is installed correctly and that the unit is at its recommended operating temperature.</li> <li>Reboot the instrument: Power OFF the instrument, wait 3 seconds, then power ON the instrument.</li> <li>If the error persists, restore the factory defaults, then re-calibrate the instrument: <ol style="list-style-type: none"> <li>Power OFF the instrument and wait 3 seconds.</li> <li>While holding down <b>Vert. Adjust</b>, power ON the instrument. This step restores the factory default settings.</li> <li>Repeat 3a–3b as needed to restore the factory defaults.</li> <li>Calibrate the vertical and horizontal axes.</li> </ol> </li> </ol>
Either of the following: <ul style="list-style-type: none"> <li>E12 or E22 errors are displayed.</li> <li>Tip Arm does not move or moves slightly.</li> </ul>	AC line voltage module is installed incorrectly.	<ol style="list-style-type: none"> <li>Determine the voltage of the electrical outlet serving the Ion OneTouch™ ES.</li> <li>Align the arrow by the correct voltage on the AC line voltage fuse module with the adjacent white arrow in the lower-right corner of the fuse socket.</li> </ol> <p>If the AC line voltage fuse module is installed incorrectly:</p> <ol style="list-style-type: none"> <li>Gently remove the module with your fingernail or a small flathead screwdriver.</li> <li>Rotate the module so that the correct voltage on the module is aligned and adjacent to the white arrow in the lower right-hand corner of the fuse socket.</li> <li>Insert the AC line voltage fuse module into the fuse socket.</li> </ol>
	Instrument is not at the recommended operating temperature	Ensure that the Ion OneTouch™ ES is at an operating temperature of 60°F to 77°F (15°C to 25°C).
Solution overflows during run	Reagent volumes are overloaded.	Repeat with reagent volumes described in enrichment procedure.
Tip is causing 8-well strip to lift out of tray slot during run	Tip is not aligned vertically.	Perform the vertical calibration procedure.
Percent template-positive ISPs after enrichment is <50% as measured by flow cytometry	Multiple causes are possible.	Contact Technical Support.

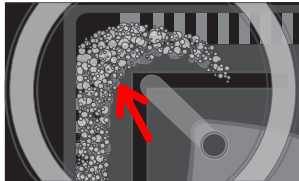
Observation	Possible cause	Recommended action
Problems with the strip position <ul style="list-style-type: none"> <li>Strip lifts up during strip push.</li> <li>Strip lifts up when tip is raised from well.</li> <li>Immediately after strip push, the strip is not in contact with the magnet.</li> </ul>	Instrument is not calibrated properly.	<ul style="list-style-type: none"> <li>Perform horizontal calibration.</li> <li>Perform vertical calibration.</li> </ul>
Tip grinds into base of instrument and Code "1999" displays	<ul style="list-style-type: none"> <li>Unit is not calibrated properly.</li> <li>Vertical calibration setting is too low or out-of-range.</li> </ul>	<ol style="list-style-type: none"> <li>Restore the factory default settings on the instrument: Hold down the vertical adjust button while powering ON the instrument. The instrument beeps several times.</li> <li>Re-calibrate the instrument.</li> <li>Perform a residual volume test.</li> </ol>
Tip hits the top of the tray at start of run	Tray is not properly seated in the instrument.	Check for debris between the tray and the instrument, then reinstall the tray. Press down firmly to ensure that tray is fully seated in the instrument.
Error messages display	Various causes are possible.	<ol style="list-style-type: none"> <li>Power the instrument OFF, then ON.</li> <li>If the error continues to display, restore the factory default settings on the instrument. Hold down the vertical adjust button while powering ON the instrument. The instrument beeps several times.</li> <li>Re-calibrate the instrument.</li> <li>Perform a residual volume test.</li> </ol>
Instrument does not aspirate or dispense liquids	Fitting(s) are loose.	<ul style="list-style-type: none"> <li>Ensure that the Luer-Lok™ connections at the elbow on the Tip Arm and at the tubing on the rear syringe pump are finger-tight.</li> <li>Ensure that the metal tip adapter fitting on the Tip Arm is finger-tight.</li> </ul> <p><b>IMPORTANT!</b> After any adjustments to the metal tip adapter, recalibrate the Ion OneTouch™ ES.</p>



## Chip Check


Observation	Possible cause	Recommended action
Chip Check fails	<ul style="list-style-type: none"> <li>Clamp is not closed.</li> <li>Chip is not properly seated.</li> <li>Debris is present on the chip socket.</li> <li>Chip is damaged.</li> </ul>	<ol style="list-style-type: none"> <li>Open the chip clamp, remove the chip, and look for signs of water outside the flow cell: <div data-bbox="971 438 1266 617" data-label="Image"> </div> </li> <li>If the chip appears damaged, replace it with a new one.</li> <li>Look for debris on the chip socket. Remove any debris by rinsing with 18-MQ water and gently dabbing the socket with a lab wipe tissue. <p><b>IMPORTANT!</b> Never rub or wipe the socket. Rubbing the socket can damage it and cause it to fail.</p> </li> <li>Close the clamp and repeat the Chip Check.</li> <li>If the chip passes, click <b>Next</b>. If the chip fails, replace it with a new chip, scan the barcode of the new chip, then press <b>Chip Check</b>.</li> <li>If Chip Check continues to fail, there could be a problem with the chip socket. Contact Technical Support.</li> </ol>

## Chip calibration (before loading sample)

Observation	Possible cause	Recommended action
Leak of unknown origin occurs	<ul style="list-style-type: none"> <li>Chip is leaking.</li> <li>Chip clamp is not closed properly.</li> <li>Problem exists with the chip clamp or socket.</li> </ul>	<ol style="list-style-type: none"> <li>Press <b>Main Menu</b>.</li> <li>Open the chip clamp, remove the chip, and gently dab the chip socket with a lab wipe tissue to absorb any fluid. <b>IMPORTANT!</b> Never rub or wipe the socket. Rubbing the socket can damage it and cause it to fail.</li> <li>Rinse the socket with 18-M<math>\Omega</math> water and gently absorb most of the water with the lab wipe.</li> <li>Repeat the rinse, then gently dab the chip socket until dry.</li> <li>Place a lab wipe on the grounding plate and dampen it with 18-M<math>\Omega</math> water. Wipe the bottom of the chip on this wipe to remove salts from the chip contacts.</li> <li>Remove the wipe, dry the grounding plate, and place chip on grounding plate. Ensure that there is no condensation outside the flow cell:</li> </ol>  <ol style="list-style-type: none"> <li>Replace the chip with a new (unused) one if needed. <b>Note:</b> The new chip can be used for sequencing after initialization completes.</li> <li>Press <b>Run</b> to restart the experiment.</li> <li>When prompted to install the new chip, ensure that the chip clamp is fully closed.</li> <li>If the chip leaks again, clean the chip socket as described above. Continued leaking, even with new chips, can indicate a chip clamp or socket problem. Contact Technical support.</li> </ol>

Observation	Possible cause	Recommended action
Error message: Calibration FAILED	<ul style="list-style-type: none"> <li>Chip is not seated in socket correctly.</li> <li>Chip is damaged.</li> </ul>	<ol style="list-style-type: none"> <li>1. Remove the chip and confirm that there is no leakage or debris on the chip socket. If leaking or debris is seen, follow the procedure for inspecting the chip and clearing debris as described under "Chip Check fails" and/or "Leak of unknown origin" above. If no leaking or debris is seen, reseal the chip in the socket.</li> <li>2. Press <b>Calibrate</b> to repeat the calibration.</li> <li>3. If the chip passes, press <b>Next</b>. If the chip still fails return to the main menu and restart the experiment with a new chip.</li> <li>4. If you continue to have chip calibration issues, there may be an issue with the chip socket. Contact Technical Support.</li> </ol>

## Chip calibration (after loading sample)

Observation	Possible cause	Recommended action
Leak of unknown origin occurs	<ul style="list-style-type: none"> <li>Chip is leaking.</li> <li>Chip clamp is not closed properly.</li> </ul>	<ol style="list-style-type: none"> <li>1. Press the <b>Abort</b> button.</li> <li>2. Open the chip clamp, remove the chip, and gently dab the chip socket with a lab wipe tissue to absorb any fluid. Do not rub or wipe the chip socket.</li> <li>3. Rinse the socket with 18 MΩ water and gently absorb most of the water with the lab wipe tissue.</li> <li>4. Repeat the rinse, then gently dab the chip socket until dry.</li> <li>5. Place a lab wipe tissue on the grounding plate and dampen it with 18 MΩ water. Wipe the bottom of the chip on this wipe to remove salts from the chip contacts.</li> <li>6. Remove the wipe, dry the grounding plate, and place the chip on the grounding plate. Check for condensation outside the flow cell:</li> </ol>  <ol style="list-style-type: none"> <li>7. If there is condensation or fluid, the chip is damaged and cannot be run.</li> <li>8. If there is no condensation or fluid, press <b>Calibrate</b> to restart the calibration procedure.</li> <li>9. If calibration passes and no leaks are visible, press <b>Next</b> to start the experiment.</li> <li>10. If the chip leaks again, clean the chip and chip socket as described above. Continued leaking can indicate a chip clamp or socket problem. Contact Technical Support.</li> </ol>




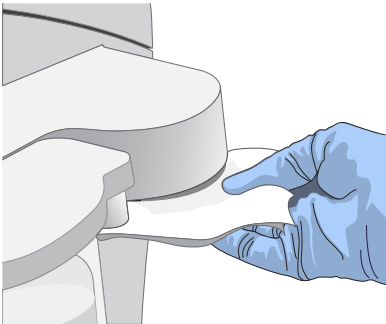
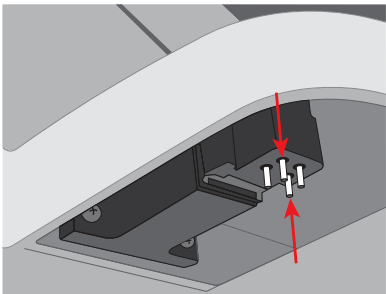
Observation	Possible cause	Recommended action
Error message: Calibration FAILED	<ul style="list-style-type: none"> <li>Chip is not seated in socket correctly.</li> <li>Chip is damaged.</li> </ul>	<ol style="list-style-type: none"> <li>1. Remove the chip and check for leaks and/or debris on the chip socket, following the procedures described in "Chip Check fails" and/or "Leak of unknown origin," above. If no leaks or debris are visible, reseal the chip in the socket.</li> <li>2. Press <b>Calibrate</b>.</li> <li>3. If the chip passes, press <b>Next</b> to start the experiment. If the chip still fails, you can try resealing the chip multiple times and pressing <b>Calibrate</b>. If you are still unable to pass calibration, press <b>Next</b> to start the run anyhow—you may still get some data on your sample.</li> <li>4. If you continue to have chip calibration issues, there may be an issue with the chip or chip socket. Contact Technical Support.</li> </ol>

## Initialization—General errors

Observation	Possible cause	Recommended action
Error message: Confirm instrument has gas pressure	Gas cylinder may be turned off or empty.	<ol style="list-style-type: none"> <li>1. Verify that the cylinder has at least 500 PSI and 30 PSI at the outlet of the regulator. Confirm that all valves between the cylinder and the Ion PGM™ Sequencer are open.</li> <li>2. Once you confirm gas pressure leading into the instrument, press <b>Yes</b> to retry verification of gas pressure. If the test continues to fail, contact Technical Support.</li> </ol>
Bottle leak check fails	<ul style="list-style-type: none"> <li>Bottle seal is not tight.</li> <li>Bottle may be damaged / defective.</li> </ul>	<ol style="list-style-type: none"> <li>1. Finger-tighten the bottles.</li> <li>2. If the bottle continues to leak, replace the bottle.</li> <li>3. If leak check continues to fail, contact Technical Support.</li> </ol>

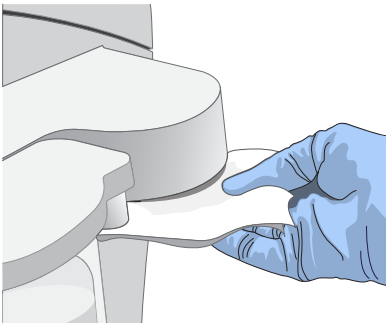
## Initialization—Auto pH errors

Observation	Possible cause	Recommended action
Error message: Please insert a chip and press <b>Start</b>	Instrument cannot detect the chip in chip socket.	<ol style="list-style-type: none"> <li>1. Open the chip clamp and remove the chip.</li> <li>2. Check for debris under the chip or in the chip socket. Remove any debris by rinsing with 18-MQ water and gently dabbing the socket with a lab wipe tissue. <b>IMPORTANT!</b> Never rub or wipe the socket. Rubbing the socket can damage it and cause it to fail.</li> <li>3. Look for liquid outside the flow cell of the chip:   </li> <li>4. If you see liquid, replace the chip with a new (unused) one. Wash the new chip once with 100% isopropanol and twice with SEQ Sample Buffer before using. <b>Note:</b> The new chip can be used for sequencing after initialization completes.</li> <li>5. Close the clamp, then press <b>Start</b> to restart the process.</li> <li>6. If the new chip also fails, there could be a problem with the chip socket. Contact Technical Support.</li> </ol>
Error message: Chip calibration failed	<ul style="list-style-type: none"> <li>• Chip is not seated in socket correctly.</li> <li>• Chip is damaged.</li> <li>• Sipper is loose.</li> </ul>	<p>Follow the procedure for "Error message: Please insert a chip and press Start."</p> <p>Follow the procedure for "Error message: Wash 2 average not stable."</p>


Observation	Possible cause	Recommended action
Error message: The system did not reach the target W2 pH and/or has a clog	The waste lines can be clogged.	<ol style="list-style-type: none"> <li>1. Press the <b>Troubleshoot</b> button. <b>Note:</b> You can skip the <b>Troubleshoot</b> button and change the chip to restart the Auto-pH routine.</li> <li>2. Remove the waste bottle.</li> <li>3. Place lab wipes under the waste arm.</li> <li>4. Gently wipe the waste arm with a lab wipe to clear liquid near the waste line.</li> </ol>  <ol style="list-style-type: none"> <li>5. Press <b>Next</b> to start buffer flow. Observe flow rates from both waste lines. One line should drip slightly faster than the other. Following the flow rate check, one of 3 results is possible: <ol style="list-style-type: none"> <li>a. If flow rate appears normal, press <b>Cancel</b> and test another chip. If Auto pH failure persists, contact Technical Support.</li> <li>b. If flow is blocked, press <b>Line Clear</b> to run the standard Line Clear procedure. If the line is unable to clear, contact Technical Support.</li> <li>c. If the result of the flow rate check is uncertain, press <b>Re-flow</b> to re-flow the buffer and re-test the flow.</li> </ol> </li> </ol> 

Observation	Possible cause	Recommended action
Error message: The system did not reach the target W2 pH ( <i>continued</i> )	Wash 1 or Wash 2 sipper may be loose.	<ol style="list-style-type: none"> <li>1. Loosen the Wash 1 cap and re-tighten the sipper. Since the gas flows when the cap is loose, tighten the sipper as quickly as possible. (The gas is not harmful to the NaOH solution and is not a hazard.)</li> <li>2. Loosen the Wash 2 cap and re-tighten the sipper. Since the gas flows when the cap is loose, tighten the sipper as quickly as possible. (The gas is not harmful to the W2 Solution and is not a hazard.)</li> <li>3. Press <b>Start</b> to re-start the auto-pH process.</li> </ol>
	Forgot to add NaOH to the Wash 1 Bottle.	<ol style="list-style-type: none"> <li>1. If there is no NaOH in the Wash 1 Bottle, loosen the cap and add 350 <math>\mu</math>L of 100 mM NaOH to the Wash 1 Bottle. (The flowing gas is not harmful to the NaOH solution and is not a hazard.)</li> <li>2. Recap the bottle and shake gently to mix.</li> <li>3. Press <b>Start</b> to restart auto-pH.</li> </ol>
	Chip is damaged.	<ol style="list-style-type: none"> <li>1. Replace the chip with a new (unused) one. Insert the chip in the socket, then press <b>Start</b>.  <b>Note:</b> The new chip can be used for sequencing after initialization completes.</li> <li>2. If the error persists, there could be a problem with the chip clamp. Contact Technical Support.</li> </ol>



Observation	Possible cause	Recommended action
Error message: W2 average not stable. Try reseating/replacing chip	Reading for W2 solution is not stabilizing quickly enough.	<ol style="list-style-type: none"> <li>1. Remove the waste bottle and gently wipe excess fluid from the waste lines with a lab wipe.</li> </ol>  <ol style="list-style-type: none"> <li>2. Check for leaks and reseat the chip (see troubleshooting for "Chip Check" and "Chip calibration" above). Replace the chip with a new (unused) one if needed. <b>Note:</b> The new chip can be used for sequencing after initialization completes.</li> <li>3. Loosen the cap in the W2 position and re-tighten the sipper. Because the gas flows when the cap is loose, tighten the sipper as quickly as possible. (The gas is not harmful and not a hazard.)</li> <li>4. After performing one or more above steps, press <b>Start</b> to re-start auto-pH. If auto-pH fails even after replacing the chip, contact Technical Support and manually adjust the pH of the Wash 2 Bottle as described in Appendix C, "Manually adjust W2 pH".</li> </ol>
Error message: W2 out of range	<ul style="list-style-type: none"> <li>• Chip measurements are very unstable.</li> <li>• Chip is damaged.</li> </ul>	See troubleshooting tips for "W2 average not stable" above.
Error message: Chip reading inconsistent. Please replace chip and try again	<ul style="list-style-type: none"> <li>• pH response of the chip is not uniform or reliable.</li> <li>• Ran out of W3 Solution or volume too low.</li> </ul>	<ol style="list-style-type: none"> <li>1. Verify that there is enough W3 Solution (&gt;25 mL) in the Wash 3 Bottle and that the sipper is secure.</li> <li>2. If necessary, loosen the Wash 3 Bottle cap, tighten the sipper, and add more W3 Solution to fill to 50 mL. Since the gas flows when the cap is loose, perform these operations as quickly as possible. (The gas is not harmful to the W3 Solution and is not a hazard.)</li> <li>3. If there is enough W3 Solution, replace the chip with a new (unused) one. Insert the chip in the socket, then press <b>Start</b>. <b>Note:</b> The new chip can be used for sequencing after initialization completes.</li> </ol>

Observation	Possible cause	Recommended action
Error message: Added too much W1 to W2	<ul style="list-style-type: none"> <li>Water quality is poor.</li> <li>18 MΩ water was exposed to air for too long.</li> <li>Incorrect solution was added to the Wash 2 Bottle.</li> <li>Too little NaOH was added to Wash 1 Bottle.</li> <li>Chip is damaged.</li> </ul>	<ol style="list-style-type: none"> <li>1. Check whether the water meets the 18 MΩ specification and 100 mM NaOH and W2 Solution were added correctly.</li> <li>2. If solutions are incorrect or water does not meet specifications, correctly prepare the solution(s) and/or use high-quality water. Abort the initialization and restart using correct solutions/water.</li> <li>3. If solutions are correct and water meets specifications, abort the initialization, return to the main menu, and proceed to the next steps.</li> <li>4. Leave the Wash 2 Bottle on the instrument.</li> <li>5. Remove the Wash 1 Bottle, leaving the sipper on the W1 port. Empty the bottle, and rinse the bottle twice with 18 MΩ water.</li> <li>6. Add 350 µL of 100 mM NaOH to the Wash 1 Bottle and reinstall on the instrument.</li> <li>7. Press <b>Initialize</b>, select the kit type, and keep pressing the <b>Next</b> button to skip all bottle prep steps until the instrument begins purging air from the bottle. Then proceed through the touchscreens as normal to complete the initialization.</li> <li>8. The next time you initialize the instrument, add 140 µL of 100 mM NaOH to the Wash 2 Bottle instead of 70 µL. Continue to use this larger volume for subsequent initializations until you receive an "Overshot Target" error message at the first auto-pH iteration, at which point follow the troubleshooting steps in "Error message: The system overshot the target W2 pH." on page 99 on the following page and then return to adding 70 µL of 100 mM NaOH.</li> <li>9. If you still receive the same initialization error ("Added too much W1 to W2"), contact Technical Support.</li> </ol>
Error message: UNDERSHOT TARGET PH: W2 pH = n.nn Failed	Auto-pH couldn't add enough Wash 1 to the Wash 2 before the maximum iterations, 10, occurred.	<ol style="list-style-type: none"> <li>1. A blockage may have occurred. Follow the procedure for "Error message: There may be a blockage or no NaOH in W1. Please check W1 and run line clear then try again."</li> <li>2. Press <b>Start</b> to re-start auto-pH. If you still get the "Undershot target pH" error, try replacing the chip with a new (unused) chip and restarting auto-pH.</li> </ol> <p><b>Note:</b> The new chip can be used for sequencing after initialization completes.</p>

Observation	Possible cause	Recommended action
Error message: The system overshot the target W2 pH.	Auto-pH added more NaOH from the Wash 1 Bottle to the Wash 2 Bottle than was needed, and reports the pH value.	<ol style="list-style-type: none"> <li>1. Press the <b>Overshoot</b> button to proceed with W2 pH adjustment.</li> <li>2. Unscrew the cap of the Wash 2 Bottle. Without removing the sipper from the bottle, lift the cap high enough to pipette 15 µL of 100 mM HCl into the Wash 2 Bottle, close and tighten cap.</li> </ol>  <ol style="list-style-type: none"> <li>3. Press <b>Next</b> to re-pressurize the Wash 2 Bottle and mix the W2 solution.</li> <li>4. Press <b>Start</b> to retry auto-pH.</li> </ol>

## Initialization—Reagent pH verification

Observation	Possible cause	Recommended action
Red failure screen, reagent pH is displayed	One or more reagents are not within the target pH.	<ol style="list-style-type: none"> <li>1. Press <b>Start</b> to repeat the pH measurements to confirm the measurement.</li> <li>2. If any reagents still fail, try replacing the chip with a new (unused) chip and repeating. <b>Note:</b> The new chip can be used for sequencing after initialization completes.</li> <li>3. If any reagents still fail, clean and re-initialize the instrument with fresh reagents and a new chip.</li> </ol>
	A possible line clog exists which persisted through the Auto pH process.	<ol style="list-style-type: none"> <li>1. From the <b>Tools</b> menu, perform a W1 line clear.</li> <li>2. Press <b>Start</b> to repeat the pH check.</li> </ol>



Observation	Possible cause	Recommended action
Red failure screen, reagent pH is <i>not</i> displayed	Chip did not calibrate.	<ol style="list-style-type: none"><li>1. Replace the chip with a new (unused) one. <b>Note:</b> The new chip can be used for sequencing after initialization completes.</li><li>2. Press <b>Start</b> to restart the pH measurement.</li><li>3. If the second test fails, contact Technical Support.</li></ol>

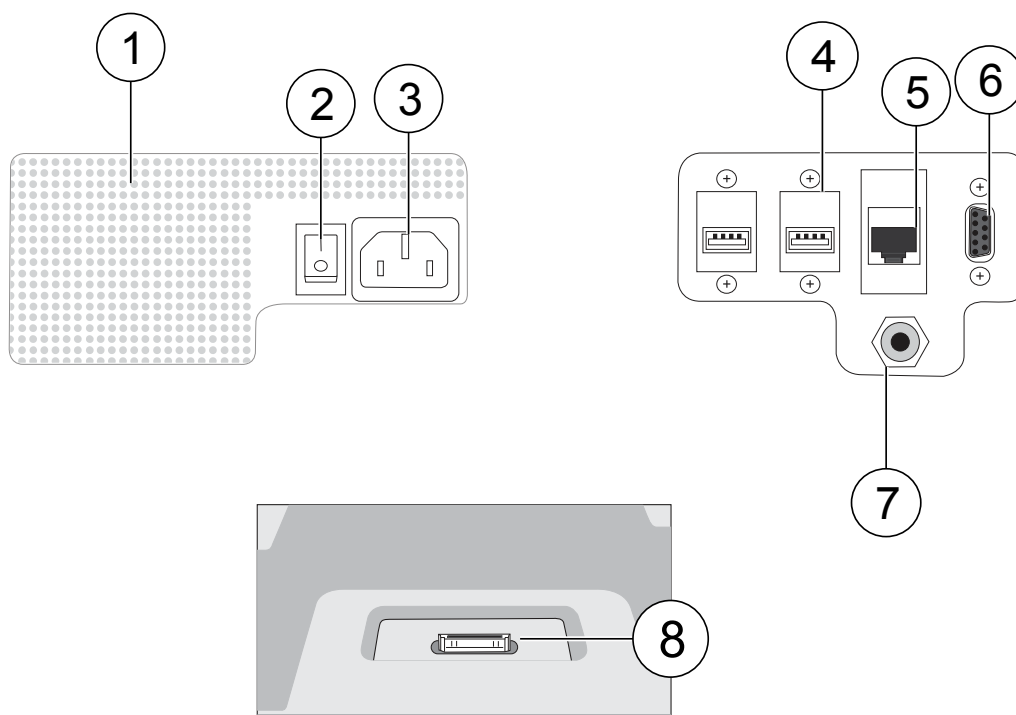
## Ion Reporter™ results

Observation	Possible cause	Recommended action
High MAPD values (>0.3) is observed.	High levels of duplicate reads occurred, which may be associated with low (<30%) polyclonality and reduced loading.	Increase library input into the isothermal amplification reaction.
	Low library read representation (<100,000 reads per sample) occurred.	Sequence poorly represented samples to a higher depth.
		Reduce the number of samples run per chip.



# Additional instrument information

## Ion PGM™ Sequencer input and output connections



Label	Component	Description
1	Instrument fan cover	<b>IMPORTANT!</b> The fan cover must be unobstructed to ensure adequate cooling and proper functioning of the Ion PGM™ Sequencer.
2	On/off switch	Power switch, where the states are on (   ) or off ( O ).
3	Power port	100-240VAC port that provides power to the instrument.
4	USB ports	Connects the barcode reader to the instrument.
5	Ethernet port	An RJ45 port that provides Ethernet (Gigabit) communication with the Ion PGM™ Sequencer.
6	RS232 port	An instrument diagnostic port

Label	Component	Description
7	Gas inlet	For nitrogen gas.
8	iPod™ port	A port for docking your iPod™ portable media player

## Power the Ion PGM™ Sequencer on or off

### Power on

**Note:** If the Ion PGM™ Sequencer is powered on, and the touchscreen is blank, touch the screen to "wake" the touchscreen.

1. Locate the power switch on the back of the instrument and turn to the on (I) position.
2. Press the power button on the front of the instrument. The switch should illuminate. When the instrument touchscreen Main Menu appears, the instrument is ready for use.
3. See "Cleaning schedule" on page 55 for when to perform 18 MΩ water or chlorite solution cleaning after powering on.

### Power off

It is not necessary to power off the instrument overnight or over the weekend. If the instrument will not be used for more than 3 days, power off the instrument as follows:

1. In the Main Menu, select **Tools > Shut Down**.
2. If you have not already cleaned the instrument, select 18 MΩ water cleaning, then press **Next** to start the cleaning process.
3. When cleaning is complete, press **Shut Down**.
4. After you exit the main touchscreen, press the **Halt** button, then **OK** when prompted. The instrument will power down.

## Update the Ion PGM™ System software

**Note:** An internet connection is required for the Ion PGM™ System to receive alerts that software updates are available.

If an update to the Ion PGM™ Sequencer software is available, the red "Alarms and Events" pop-up appears in the touchscreen Main Menu to alert you. Press the red pop-up to see the detailed messages. If a message states New Software Available, update the software as follows:

1. In the Main Menu, select **Options ▶ Updates**.
2. Select the **Released Updates** checkbox, then press **Check**.

3. When the message Press Update to begin update process appears, press **Update**.

**Note:** If the message All Software Current appears, press **Back** to return to the Main Menu.

4. When the message Installing Completed displays, follow the onscreen prompts to restart the instrument.

**Note:** In some cases, the instrument restarts automatically after software installation.



# Manually adjust W2 pH

## Materials and equipment needed

- Orion™ 3-Star Plus pH Benchtop Meter Kit or equivalent
- Nitrogen gas tank, tube, and flow meter
- 100 mM NaOH (prepared fresh daily)
- Pipette tips and pipette
- Magnetic stirrer and stir bar
- 100 mM HCl

## Procedure

If an error message during the automatic pH process indicates that there is a problem adjusting the pH of the W2 Solution, use the following procedure to adjust the pH of the W2 Solution in the Wash 2 Bottle manually.

1. Before proceeding, rinse an empty Wash 2 Bottle and have it ready next to the instrument. Also have an extra Wash 2 Bottle cap ready.  
**Note:** Gas will be flowing out of the Wash 2 cap, so perform the next steps as quickly as possible (flowing gas will not harm the W2 Solution, and is not a hazard).
2. Remove the Wash 2 Bottle attached to the instrument, then cap the bottle.
3. Secure the empty Wash 2 Bottle (from step 1) to the instrument—do not remove the sipper. This bottle contains the gas flowing out of the instrument while adjust the pH of the W2 Solution, and protects the sipper from contamination.
4. Move the Wash 2 Bottle containing the W2 Solution to the stir plate near the nitrogen gas tube.
5. Secure the gas tube so that it extends inside the mouth of the Wash 2 Bottle but not below the surface of the W2 Solution.
6. Set the gas flow to 0.5 lpm. Start mixing the W2 Solution fast sufficient for a small whirlpool to form.
7. Calibrate the pH meter using a three-point calibration. Rinse any buffering solution from the pH probe before preparing solutions.





8. Adjust the pH of the W2 Solution to  $7.55 \pm 0.1$  by adding a small amount of freshly prepared 100 mM NaOH to the solution, then measuring the pH using the pH meter. Add small aliquots, then allow the pH to equilibrate before adding more.  
**Note:** If the pH rises above 7.75, use 100 mM hydrochloric acid (HCl) to readjust the pH to  $7.55 \pm 0.1$ .
9. When the pH is stable, turn off the gas, remove the gas line, then cap the Wash 2 Bottle.
10. Move the bottle to the instrument, remove the empty Wash 2 Bottle from the instrument, then place the sipper inside the Wash 2 Bottle whose pH adjusted.
11. Secure the cap firmly. Press **Next** to exit the automated pH check, then continue with instrument initialization.



# Safety



**WARNING! GENERAL SAFETY.** Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.




- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
- Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, etc). To obtain SDSs, see the “Documentation and Support” section in this document.

## Symbols on this instrument

Symbols may be found on the instrument to warn against potential hazards or convey important safety information. In this document, the hazard symbol is used along with one of the following user attention words:



- **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, will result in death or serious injury.

Symbol	English	Français
	Caution, risk of danger Consult the manual for further safety information.	Attention, risque de danger Consulter le manuel pour d'autres renseignements de sécurité.
	Protective conductor terminal (main ground)	Borne de conducteur de protection (mise à la terre principale)

Symbol	English	Français
	<p>Do not dispose of this product in unsorted municipal waste</p> <p> <b>CAUTION!</b> To minimize negative environmental impact from disposal of electronic waste, do not dispose of electronic waste in unsorted municipal waste. Follow local municipal waste ordinances for proper disposal provision and contact customer service for information about responsible disposal options.</p>	<p>Ne pas éliminer ce produit avec les déchets usuels non soumis au tri sélectif.</p> <p> <b>CAUTION!</b> Pour minimiser les conséquences négatives sur l'environnement à la suite de l'élimination de déchets électroniques, ne pas éliminer ce déchet électronique avec les déchets usuels non soumis au tri sélectif. Se conformer aux ordonnances locales sur les déchets municipaux pour les dispositions d'élimination et communiquer avec le service à la clientèle pour des renseignements sur les options d'élimination responsable.</p>

## Safety alerts on this instrument

Additional text may be used with one of the symbols described above when more specific information is needed to avoid exposure to a hazard. See the following table for safety alerts found on the instrument.

English	French translation
<p> <b>CAUTION! Hazardous chemicals.</b> Read the Safety Data Sheets (SDSs) before handling.</p>	<p><b>ATTENTION! Produits chimiques dangereux.</b> Lire les fiches signalétiques (FS) avant de manipuler les produits.</p>
<p> <b>CAUTION! Hazardous waste.</b> Refer to SDS(s) and local regulations for handling and disposal.</p>	<p><b>ATTENTION! Déchets dangereux.</b> Lire les fiches signalétiques (FS) et la réglementation locale associées à la manipulation et à l'élimination des déchets.</p>



## Safety information for instruments not manufactured by Thermo Fisher Scientific

Some of the accessories provided as part of the instrument system are not designed or built by Thermo Fisher Scientific. Consult the manufacturer's documentation for the information needed for the safe use of these products.

### Instrument safety

#### General



**CAUTION! Do not remove instrument protective covers.** If you remove the protective instrument panels or disable interlock devices, you may be exposed to serious hazards including, but not limited to, severe electrical shock, laser exposure, crushing, or chemical exposure.

#### Physical injury



**CAUTION! Moving Parts.** Moving parts can crush, pinch and cut. Keep hands clear of moving parts while operating the instrument. Disconnect power before servicing.

#### Electrical



**WARNING! Ensure appropriate electrical supply.** For safe operation of the instrument:

- Plug the system into a properly grounded receptacle with adequate current capacity.
- Ensure the electrical supply is of suitable voltage.
- Never operate the instrument with the ground disconnected. Grounding continuity is required for safe operation of the instrument.



**WARNING! Power Supply Line Cords.** Use properly configured and approved line cords for the power supply in your facility.



**WARNING! Disconnecting Power.** To fully disconnect power either detach or unplug the power cord, positioning the instrument such that the power cord is accessible.

## Cleaning and decontamination



**CAUTION! Cleaning and Decontamination.** Use only the cleaning and decontamination methods specified in the manufacturer's user documentation. It is the responsibility of the operator (or other responsible person) to ensure the following requirements are met:

- No decontamination or cleaning agents are used that could cause a HAZARD as a result of a reaction with parts of the equipment or with material contained in the equipment.
- The instrument is properly decontaminated a) if hazardous material is spilled onto or into the equipment, and/or b) prior to having the instrument serviced at your facility or sending the instrument for repair, maintenance, trade-in, disposal, or termination of a loan (decontamination forms may be requested from customer service).
- Before using any cleaning or decontamination methods (except those recommended by the manufacturer), users should confirm with the manufacturer that the proposed method will not damage the equipment.

## Laser



**CAUTION! LASER HAZARD, Bar Code Scanner.** The bar code scanner included with the instrument system is a Class 2 laser. To avoid damage to eyes, do not stare directly into the beam or point into another person's eyes.

## Safety and electromagnetic compatibility (EMC) standards

The instrument design and manufacture complies with the standards and requirements for safety and electromagnetic compatibility as noted in the following table:

### Safety

Reference	Description
EU Directive 2006/95/EC	European Union "Low Voltage Directive"
IEC 61010-1 EN 61010-1 UL 61010-1 CSA C22.2 No. 61010-1	<i>Safety requirements for electrical equipment for measurement, control, and laboratory use – Part 1: General requirements</i>
IEC 61010-2-010 EN 61010-2-010	<i>Safety requirements for electrical equipment for measurement, control and laboratory use – Part 2-010: Particular requirements for laboratory equipment for the heating of materials</i>

## EMC

Reference	Description
Directive 2004/108/EC	European Union “EMC Directive”
EN 61326-1	<i>Electrical Equipment for Measurement, Control and Laboratory Use – EMC Requirements – Part 1: General Requirements</i>
FCC Part 15	U.S. Standard “Industrial, Scientific, and Medical Equipment”
AS/NZS 2064	<i>Limits and Methods of Measurement of Electromagnetic Disturbance Characteristics of Industrial, Scientific, and Medical (ISM) Radiofrequency Equipment</i>
ICES-001, Issue 3	<i>Industrial, Scientific and Medical (ISM) Radio Frequency Generators</i>

## Environmental design

Reference	Description
Directive 2012/19/EU	European Union “WEEE Directive” – Waste electrical and electronic equipment
Directive 2011/65/EU	European Union “RoHS Directive” – Restriction of hazardous substances in electrical and electronic equipment

## Chemical safety



**WARNING! GENERAL CHEMICAL HANDLING.** To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below. Consult the relevant SDS for specific precautions and instructions:

- 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. To obtain SDSs, see the “Documentation and Support” section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of 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.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the 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



**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. Conduct all work in properly equipped facilities with the appropriate safety equipment (for example, physical containment devices). Safety equipment can also include items for personal protection, such as gloves, coats, gowns, shoe covers, boots, respirators, face shields, safety glasses, or goggles. Individuals should be trained according to applicable regulatory and company/ institution requirements before working with potentially biohazardous materials. Follow all applicable local, state/provincial, and/or national regulations. The following references provide general guidelines when handling biological samples in laboratory environment.

- U.S. Department of Health and Human Services, *Biosafety in Microbiological and Biomedical Laboratories (BMBL)*, 5th Edition, HHS Publication No. (CDC) 21-1112, Revised December 2009; found at:  
[www.cdc.gov/biosafety/publications/bmbl5/BMBL.pdf](http://www.cdc.gov/biosafety/publications/bmbl5/BMBL.pdf)
- World Health Organization, *Laboratory Biosafety Manual*, 3rd Edition, WHO/CDS/CSR/LYO/2004.11; found at:  
[www.who.int/csr/resources/publications/biosafety/Biosafety7.pdf](http://www.who.int/csr/resources/publications/biosafety/Biosafety7.pdf)



# Documentation and support

## Customer and technical support

Visit **[thermofisher.com/support](http://thermofisher.com/support)** for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
  - Product FAQs
  - Software, patches, and updates
  - Training for many applications and instruments
- Order and web support
- Product documentation, including:
  - User guides, manuals, and protocols
  - Certificates of Analysis
  - Safety Data Sheets (SDSs; also known as MSDSs)

**Note:** For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at **[www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html)**. If you have any questions, please contact Life Technologies at **[www.thermofisher.com/support](http://www.thermofisher.com/support)**.

For support visit [thermofisher.com/support](http://thermofisher.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)  
[thermofisher.com](http://thermofisher.com)

7 November 2016

**ThermoFisher**  
S C I E N T I F I C