

SuperScript™ VILO™ MasterMix

Catalog Number 11755-050, 11755-250, and 11755-500

Doc. Part No. 100012386 Pub. No. MAN0004286 Rev. A.0



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product description

Invitrogen™ SuperScript™ VILO™ MasterMix provides the high-temperature capability of SuperScript™ III Reverse Transcriptase (RT) in an optimized format for generating first-strand cDNA for use in real-time quantitative RT-PCR (qRT-PCR). This formulation can be used with very low and very high amounts of input RNA (up to 2.5 µg total RNA in a 20-µL reaction).

SuperScript™ VILO™ Master Mix includes SuperScript™ III RT, RNaseOUT™ Recombinant Ribonuclease Inhibitor, a proprietary helper protein, random primers, MgCl₂, and dNTPs.

SuperScript™ III RT is an engineered version of M-MLV RT with reduced RNase H activity and increased thermal stability. The enzyme can be used to synthesize cDNA at a temperature range of 42–55°C. Because SuperScript™ III RT is not significantly inhibited by ribosomal and transfer RNA, it can be used to synthesize cDNA from total RNA. RNaseOUT™ Recombinant Ribonuclease Inhibitor safeguards against the degradation of target RNA due to ribonuclease contamination.

Contents and storage

Contents	Cat. no. 11755-050 (50 reactions)	Cat. no. 11755-250 (250 reactions)	Cat. no. 11755-500 (500 reactions)	Storage conditions
SuperScript™ VILO™ Master Mix	200 µL	1000 µL	2000 µL	Store at -20°C (non-frost-free)

Important guidelines

cDNA synthesis guidelines

- High-quality, intact RNA is essential for accurate quantification in qPCR. RNA should be devoid of RNase contamination and aseptic conditions should be maintained. RNA quality can be analyzed using a bioanalyzer or by agarose gel electrophoresis.
- Starting material can range up to 2.5 µg of total RNA in a 20-µL cDNA synthesis reaction. RNA quantity can be determined using UV absorbance at 260 nm or the Qubit™ RNA Assay Kit and Qubit™ 2.0 Fluorometer (see “Related products” on page 4 for ordering information).
- To isolate total RNA, we recommend TRIzol™ Reagent, the PureLink™ RNA Mini Kit, or the MagMAX™-96 Total RNA Isolation Kit (see “Related products” on page 4). Isolation of mRNA is typically not necessary, although incorporating this step may improve the yield of specific cDNAs.
- DNase I, Amplification Grade, may be used to eliminate genomic DNA contamination from the total RNA (see “Related products” on page 4).
- Shorter incubation times and/or higher temperatures may be used (e.g., 50°C for 30 minutes), but may result in reduced yields of cDNA.
- For increased yields of cDNA, longer incubation times may be used (up to 120 minutes at 42°C).

qPCR using fluorescent primers or probes

Up to 10% of the qPCR reaction volume may be undiluted cDNA (e.g., for a 20-µL qPCR, use up to 2 µL of undiluted cDNA).

qPCR using SYBR™ Green or SYBR™ GreenER™ reagent

If you started with ≤100 ng of total RNA, up to 10% of the qPCR reaction volume may be undiluted cDNA (e.g., for a 20-µL qPCR, use up to 2 µL of undiluted cDNA).

If you started with >100 ng total RNA, we recommend testing a serial dilution of cDNA in qPCR for optimal results. Higher concentrations of cDNA may affect the signal baseline in SYBR™ Green and SYBR™ GreenER™ reactions.

Methods

Synthesize first-strand cDNA

The following protocol has been optimized for generating first-strand cDNA for use in two-step qRT-PCR. The reaction volume may be scaled as needed up to 100 μL . A negative RT control protocol is provided below.

1. For a single reaction, combine the following components in a sterile PCR tube or plate well on ice.

Component	Volume
SuperScript™ VILO™ MasterMix	4 μL
RNA (up to 2.5 μg)	X μL
DEPC-treated water	to 20 μL

2. Gently mix and incubate at 25°C for 10 minutes.
3. Incubate at 42°C for 60 minutes.
4. Terminate the reaction at 85°C at 5 minutes.
5. Use the diluted or undiluted cDNA in qPCR or store at -20°C.

Prepare negative RT control

1. For a volume of RNA = X μL , add the following to a sterile PCR tube or plate well on ice.

Component	Volume
SuperScript™ VILO™ MasterMix	4 μL
DEPC-treated water	16 - X μL

2. Incubate at 65°C for 10 minutes to denature the reverse transcriptase.
3. Add X μL of RNA (up to 2.5 μg) for a total reaction volume of 20 μL .
4. Proceed with steps 2-5 from "Synthesize first-strand cDNA" on page 3.

Related products

Product	Amount	Cat. no.
TRIzol™ Reagent	100 mL	15596-026
	200 mL	15596-018
PureLink™ RNA Mini Kit	10 preps	12183020
	50 preps	12183018A
MagMAX™-96 Total RNA Isolation Kit	96 rxns	AM1830
Qubit™ RNA Assay Kit	500 assays	Q32855
Qubit™ 2.0 Fluorometer	1 unit	Q32866
DNase I, Amplification Grade	100 units	18068-015

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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

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.

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 | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2015 Thermo Fisher Scientific Inc. All rights reserved. TRIzol is a trademark of Molecular Research Center, Inc.

For support visit thermofisher.com/support or email techsupport@lifetech.com

thermofisher.com

2 November 2015

ThermoFisher
SCIENTIFIC