

M-MLV Reverse Transcriptase

Catalog Number 28025-013 and 28025-021

Doc. Part No. 28025.PPS Pub. No. MAN0001462 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

Moloney Murine Leukemia Virus Reverse Transcriptase (M-MLV RT) uses single-stranded RNA or DNA in the presence of a primer to synthesize a complementary DNA strand. This enzyme is isolated from *E. coli* expressing a portion of the *pol* gene of M-MLV on a plasmid. The enzyme is used to synthesize first-strand cDNA up to 7 kb.

Contents and storage

Component	Quantity	Storage conditions
M-MLV RT (200 U/ μ L)	40,000 units (Cat. no. 28025-013) 200,000 units (Cat. no. 28025-021)	<ul style="list-style-type: none"> Store at -20°C (non-frost-free). Refreeze 5X First-Strand Buffer and 0.1 M DTT immediately after use.
5X First-Strand Buffer ^[1]	1 mL	
DTT	0.1 M	

^[1] 250 mM Tris-HCl (pH 8.3 at room temperature), 375 mM KCl, 15 mM MgCl₂

Required materials not provided

RNaseOUT™ Recombinant Ribonuclease Inhibitor (40 units/ μ L; Cat. no. 10777-019).

Before you begin

- Prepare storage buffer: 20 mM Tris-HCl (pH 7.5), 100 mM NaCl, 0.1 mM EDTA, 1 mM DTT, 0.01% (v/v) NP-40, 50% (v/v) glycerol
- Thaw 5X First-Strand Buffer and 0.1 M DTT at room temperature just before use
Note: Refreeze immediately after use.

For Research Use Only. Not for use in diagnostic procedures.

Synthesize first-strand cDNA using M-MLV RT

A 20- μL reaction volume can be used for 1 ng–5 μg of total RNA or 1–500 ng of mRNA.

1. Add the following components to a nuclease-free microcentrifuge tube:

Oligo (dT) ₁₂₋₁₈ (500 $\mu\text{g}/\text{mL}$), or 50–250 ng random primers, or 2 pmole gene-specific primer	1 μL
Total RNA, or mRNA	1 ng to 5 μg total RNA, or 1 ng to 500 ng of mRNA
10 mM dNTP Mix (10 mM each dATP, dGTP, dCTP and dTTP at neutral pH)	1 μL
Sterile, distilled water	To 12 μL

2. Heat mixture to 65°C for 5 minutes and quick chill on ice. Collect the contents of the tube by brief centrifugation and add:

5X First-Strand Buffer	4 μL
0.1 M DTT	2 μL
RNaseOUT™ Recombinant Ribonuclease Inhibitor (40 units/ μL)	1 μL

Note: When using less than 50 ng of starting RNA, the addition of RNaseOUT™ is essential.

3. Mix contents of the tube gently and incubate at 37°C for 2 minutes.
4. Add 1 μL (200 units) of M-MLV RT, and mix by pipetting gently up and down. If using random primers, incubate tube at 25°C for 10 minutes.

Note: If less than 1 ng of RNA is used, reduce the amount of M-MLV RT in the reaction to 0.25 μL (50 units), and add the sterile, distilled water to 20 μL final volume.

5. Incubate 50 minutes at 37°C.
6. Inactivate the reaction by heating at 70°C for 15 minutes.

The cDNA can now be used as a template for amplification in PCR. However, amplification of some PCR targets (>1 kb) may require the removal of RNA complementary to the cDNA. To remove RNA complementary to the cDNA, add 1 μL (2 units) of *E. coli* RNase H and incubate at 37°C for 20 minutes.

Prepare PCR reaction

Use only 10% of the first-strand reaction (2 μL of the reaction from “Synthesize first-strand cDNA using M-MLV RT” on page 2) for PCR. Adding larger amounts of the first-strand reaction may not increase amplification and may result in decreased amounts of PCR product.

1. Add the following to a PCR reaction tube for a final reaction volume of 50 μL :

10X PCR Buffer [200 mM Tris-HCl (pH 8.4), 500 mM KCl]	5 μL
50 mM MgCl_2 ^[1]	1.5 μL
Amplification primer 1 (10 μM)	1 μL
Amplification primer 2 (10 μM)	1 μL
Taq DNA polymerase (5 U/ μL)	0.4 μL
cDNA ^[2]	2 μL
Autoclaved, distilled water	38.1 μL

^[1] For best results, determine the optimal concentration of MgCl_2 empirically for each template-primer pair.

^[2] From “Synthesize first-strand cDNA using M-MLV RT” on page 2.

2. Mix gently and layer 1–2 drops (~50 μL) of silicone oil over the reaction.

Note: The addition of silicone oil is unnecessary in thermal cyclers equipped with a heated lid.

3. Heat reaction to 94°C for 2 minutes to denature.
4. Perform 15 to 40 cycles of PCR. Annealing and extension conditions are primer and template dependent and must be determined empirically.

Unit definition

One unit incorporates 1 nmole of dTTP into acid-precipitable material in 10 minutes at 37°C using poly(A)•oligo(dT)₂₅ as template-primer.

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

©2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

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

thermofisher.com

30 October 2015

ThermoFisher
SCIENTIFIC