

Maxima Reverse Transcriptase

Catalog Number EP0741, EP0742, EP0743

Pub. No. MAN0012044 Rev. D.00



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.

Contents and storage

Cat No.	Contents	Amount	Storage
EP0741	Maxima Reverse Transcriptase, 200 U/μL	2000 U	-25 °C to -15 °C
	5X RT Buffer*	1 mL	
EP0742	Maxima Reverse Transcriptase, 200 U/μL	10000 U	
	5X RT Buffer*	2 x 1 mL	
EP0743	Maxima Reverse Transcriptase, 200 U/μL	4 x 10000 U	
	5X RT Buffer*	4 x 1 mL	

*5X RT Buffer is also available separately (#B91)

Description

Thermo Scientific™ Maxima™ Reverse Transcriptase (RT) is a novel reverse transcription enzyme that was developed by Thermo Scientific through *in vitro* evolution of M-MuLV RT. The enzyme possesses an RNA and DNA-dependent polymerase activity as well as RNase H activity.

Features

- High yields of full-length cDNA up to 20 kb
- Active up to 65 °C
- Thermostable – 90% active after incubation at 50 °C for 60 minutes in a reaction mixture
- High sensitivity - reproducible cDNA synthesis from a wide range of starting total RNA amounts (1 pg - 5 μg)
- Efficient – complete cDNA synthesis in 15-30 minutes
- Incorporates modified nucleotides

Applications

- First strand cDNA synthesis
- RT-PCR
- RT-qPCR
- DNA labeling
- Primer extension

Source

E. coli cells carrying an engineered *pol* gene fragment of Moloney Murine Leukemia Virus.

Definition of Activity Unit

One unit of the enzyme incorporates 1 nmol of dTMP into a polynucleotide fraction in 10 min at 37 °C.

Storage Buffer

The enzyme is supplied in: 50 mM Tris-HCl (pH 7.5), 0.1 M NaCl, 1 mM EDTA, 5 mM DTT, detergent and 50% (v/v) glycerol.

5X RT Buffer

250 mM Tris-HCl (pH 8.3 at 25 °C), 375 mM KCl, 15 mM MgCl₂, 50 mM DTT.

Inhibition and Inactivation

- Inhibitors: metal chelators, inorganic phosphate, pyrophosphate, and polyamines.
- Inactivated by heating at 85 °C for 5 min.

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

Protocol for First Strand cDNA Synthesis

The following protocol is optimized to generate first-strand cDNA for use in two-step RT-PCR.

Mix and briefly centrifuge all components after thawing, keep on ice.

1. Add reaction components into a sterile, nuclease-free tube on ice in the indicated order:

Components	Volume
Template RNA: total RNA poly(A) RNA specific RNA	1 pg - 5 µg 0.1 pg - 500 ng 0.01 pg - 500 ng
Primer: Oligo(dT) ₁₈ (#SO131) Random hexamer (#SO142) gene-specific primer	1 µL (100 pmol) 1 µL (100 pmol) 15-20 pmol
dNTP Mix, 10 mM each (#R0191)	1 µL (0.5 mM final concentration)
Water, nuclease-free	to 14.5 µL

2. **Optional:** If the RNA template is GC-rich or is known to contain secondary structures, mix gently, centrifuge briefly and incubate at 65 °C for 5 min. Chill on ice, briefly centrifuge again and place on ice.

3. Add the following components in the indicated order:

Components	Volume
5X RT Buffer	4 µL
Thermo Scientific™ RiboLock RNase Inhibitor (#EO0381)	0.5 µL (20 U)
Maxima Reverse Transcriptase	1 µL (200 U)
Total volume	20 µL

Mix gently and centrifuge briefly.

4. Incubate:

- if an oligo(dT)₁₈ primer or gene-specific primer is used, incubate for 30 min at 50 °C.
- if a random hexamer primer is used, incubate for 10 min at 25 °C followed by 30 min at 50 °C.
For transcription of GC-rich RNA, the reaction temperature can be increased to 65 °C.

5. Terminate the reaction by heating at 85 °C for 5 minutes.

Note

- The reverse transcription reaction product can be used directly in PCR or qPCR, or stored at -20 °C for up to one week. For longer storage, -70 °C is recommended. Avoid freeze/thaw cycles of the cDNA.
- Use 2 µL of the reaction mix to perform PCR in a 50 µL volume.

Recommendations for two-step RT-qPCR

- **Priming:** use a mix of oligo (dT)₁₈ and random primers 25 pmol each per 20 µL reaction.
- **Incubation:** 10 min at 25 °C followed by 15 min at 50 °C.

Recommendations for long RT-PCR (>5 kb)

- **Priming:** oligo (dT)₁₈ or gene specific primer should be used.
- Use 20 U of Maxima Reverse Transcriptase per reaction. 1X RT buffer can be used to dilute the enzyme just prior to reaction.
- **Incubation:** 30 min at 50 °C.

Revision history: Pub. No. MAN0012044

Revision	Date	Description
D.00	2024-01-18	Revized user guide template, removed COA content and updated storage buffer composition.

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 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.



Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 Vilnius, Lithuania
For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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 this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

18 January 2024

ThermoFisher
SCIENTIFIC