

Maxima H Minus cDNA Synthesis Master Mix

Catalog Number M1661, M1662

Pub. No. MAN0016392 Rev. C00



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](https://www.thermofisher.com/support).

Contents

Cat. No.	Contents	Amount	Storage
M1661 (50 rxns)	Maxima H Minus cDNA Synthesis Master Mix (5X)	200 μ L	-25 °C to -15 °C
	Maxima No RT Control	200 μ L	
	Water, nuclease-free	1.25 mL	
M1662 (200 rxns)	Maxima H Minus cDNA Synthesis Master Mix (5X)	800 μ L	
	Maxima No RT Control	200 μ L	
	Water, nuclease-free	1.25 mL	

Description

The Thermo Scientific™ Maxima™ H Minus cDNA Synthesis Master Mix provides a simple workflow that allows cDNA synthesis in a one-tube procedure. The system is convenient, and optimized for cDNA synthesis in two-step real time quantitative RT-PCR (RT-qPCR) applications.

The master mix contains Maxima H Minus Reverse Transcriptase (RT) and Thermo Scientific™ RiboLock™ RNase Inhibitor. Maxima H Minus Reverse Transcriptase is an advanced enzyme derived from M-MuLV RT. The enzyme features high thermostability, robustness, and increased cDNA synthesis rate compared to wild type M-MuLV RT. The recombinant RiboLock RNase Inhibitor effectively protects RNA template from degradation by RNases A, B, and C at temperatures up to 55 °C. The master mix also contains reaction buffer, dNTPs, oligo (dT)₁₈, and random hexamer primers.

The Maxima H Minus cDNA Synthesis Master Mix is capable of reproducible cDNA synthesis starting with 1 pg to 5 μ g of total RNA at elevated temperatures (50–65 °C). The synthesis reaction can be completed in 15–30 minutes.

PROCEDURAL GUIDELINES

Avoiding ribonuclease contamination

RNA purity and integrity is essential for synthesis of full-length cDNA. RNA can be degraded by RNase A, which is a highly stable contaminant found in laboratory environments. General recommendations to avoid RNase contamination:

- Use certified nuclease-free labware or DEPC-treat all tubes and pipette tips to be used in cDNA synthesis.
- Wear gloves when handling RNA and reagents, as skin is a common source of RNases. Change gloves frequently.
- Use RNase-free reagents, including high quality water (e.g., Water, nuclease-free, #R0581).
- Keep the kit components tightly sealed when not in use. Keep all tubes tightly closed during the reverse transcription reaction.

Template RNA

- The kit is compatible with total cellular RNA isolated by standard methods. Purified RNA must be free of salts, metal ions, ethanol, and phenol to avoid inhibiting the cDNA synthesis reaction.
- For RT-qPCR applications, template RNA must be free of DNA contamination. Prior to cDNA synthesis, treat RNA samples with DNase I, RNase-free ([#EN0521](#)) or dsDNase ([#EN0771](#)) to remove residual DNA (See product pages at www.thermofisher.com for detailed protocols).
- Always perform a No RT Control reaction, which includes all components for reverse transcription except the Maxima H Minus RT.

RNA sample quality

- Assess RNA integrity prior to cDNA synthesis. Total eukaryotic RNA can be analyzed by agarose gel electrophoresis followed by ethidium bromide staining.
- The RNA is considered intact, if both 18S and 28S rRNA appear as sharp bands after electrophoresis of total RNA. The 28S rRNA band should be approximately twice as intense as the 18S rRNA. Smearing of rRNA bands is an indication of degraded mRNA, and a new sample of total RNA should be prepared.
- Alternatively, total RNA can be analyzed using a bioanalyzer (e.g., Agilent 2100) to obtain a RNA integrity number, which provides quantitative information about the general state of the RNA sample (2). A reference gene/target gene 3':5' integrity assay (3) can also be used to determine the integrity of the RNA sample.

Primers

- The Maxima H Minus cDNA Synthesis Master Mix contains oligo(dT)₁₈ and random hexamer primers to prime synthesis of first strand cDNA. This primer mixture ensures high sensitivity in low copy number transcript detection assays.

PROTOCOL

Allow kit components to thaw, then mix and briefly centrifuge. Store the tubes on ice.

1. Add the following reagents into a sterile, RNase free tube on ice in the following order:

Component	Volume
Maxima H Minus cDNA Synthesis Master Mix (5X)	4 µL
Template RNA Total RNA or poly(A) mRNA or Specific RNA	1 pg to 5 µg 0.1 pg to 500 ng 0.01 pg to 500 ng
Water, nuclease-free	to 20 µL
Total volume	20 µL

2. Mix gently and centrifuge.
3. Incubate at 25 °C for 10 minutes.
4. Incubate at 50 °C for 15 minutes. **Note.** If using >1 µg RNA template, increase the reaction time to 30 minutes. For RNA templates that are GC-rich or have a large amount of secondary structure, the reaction temperature can be increased to 65 °C.
5. Terminate the reaction by heating at 85 °C for 5 minutes.

The reaction product of the first strand cDNA synthesis can be used directly in qPCR. Store at –20 °C for up to one week, or –70 °C for long term storage. Avoid freeze-thaw cycles of the cDNA.

qPCR

Use the product of the cDNA synthesis reaction directly in qPCR. Normally, 2 µL of the RT product is used as template for subsequent qPCR in a 25 µL total volume.

CONTROL REACTIONS

Use the following negative control reactions to verify the results of the first strand cDNA synthesis.

- **No RT control** to assess genomic DNA contamination of the RNA sample. The Maxima No RT Control reaction includes all reagents for the reverse transcription reaction except the Maxima H Minus RT.
- **No template control (NTC)** to assess for reagent contamination. The NTC reaction includes all reagents necessary for the reverse transcription reaction except the RNA template.

Prepare control reactions according to the following instructions.

Note: Template RNA amount and reverse transcription incubation conditions should reflect the conditions used for experimental samples.

1. Add the following reagents into a sterile, RNase free tube on ice in the following order:

Component	No RT control
Maxima H Minus cDNA Synthesis Master Mix (5X)	—
Template RNA	varies
Water, nuclease-free	to 20 µL
Total volume	20 µL

2. Mix gently and centrifuge.

3. Incubate at 25 °C for 10 minutes.

4. Incubate at 50 °C for 15 minutes.

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

Troubleshooting

No qPCR product generated or product appears late in qPCR
<p>Poor integrity of RNA template RNA purity and integrity is essential for synthesis and quantification of cDNA.</p> <ul style="list-style-type: none"> • Always assess the integrity of RNA prior to cDNA synthesis. • Use freshly prepared RNA. • Multiple freeze/thaw cycles of the RNA sample and synthesized cDNA is not recommended. • Follow general recommendations to avoid RNase contamination and discard low quality RNA. <p>Low template purity (inhibitors in RNA sample) Trace amounts of agents used in RNA purification protocols may remain in solution and inhibit first strand synthesis (e.g., SDS, EDTA, guanidine salts, phosphate, pyrophosphate, polyamines, spermidine). Remove trace contaminants by re-precipitating the RNA with ethanol and wash the pellet with 75 % ethanol.</p> <p>Insufficient template quantity Increase the amount of RNA template in the first strand reaction to the recommended level.</p> <p>GC-rich template If the RNA template is GC-rich or is known to contain secondary structures, the temperature of the reverse transcription reaction can be increased up to 65 °C.</p> <p>Excess amount of cDNA in qPCR Decrease amount of cDNA synthesis reaction in qPCR. The volume of the cDNA synthesis reaction mixture should not exceed 10 % of the final PCR reaction mixture.</p>
RT-qPCR product in No RT Control
<p>RNA template is contaminated with DNA. The presence of a PCR product in the negative control (No RT Control) reaction indicates that the reaction is contaminated with DNA. Follow the protocol carefully and make sure that dsDNase treatment step prior reverse transcription is not omitted. To further enhance genomic DNA elimination efficiency, template RNA incubation with dsDNase step can be prolonged to 5 minutes.</p>

References

1. Wiame, I., et al., Irreversible heat inactivation of DNaseI without RNA degradation, *BioTechniques*, 29, 252-256, 2000.
2. Fleige, S., Pfaffl, M.W., RNA integrity and the effect on the real-time qRT-PCR performance, *Mol. Aspects Med.*, 27, 126-139, 2006.
3. Nolan, T., et al., Quantification of mRNA using real-time RT-PCR, *Nat. Protoc.*, 1, 1559-1582, 2006.

Revision history: Pub. No. MAN0016392

Revision	Date	Description
C00	2024-02-26	Revized user guide template

Limited product warranty

Life Technologies Corporation and/or it 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.