


High-Capacity RNA-to-cDNA™ Kit

Catalog Numbers 4387406 (50 reactions), 4388950 (500 reactions)

Pub. No. 4387949 Rev. D

 **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

Upon receipt, store the kit at -25°C to -15°C for maximum shelf life.

Contents	50 reactions	500 reactions	Storage
2X RT Buffer Mix (50 × 20-µL reactions)	500 µL	5 × 1000 µL	Store at 2°C to 8°C for up to 6 months, or -25°C to -15°C for long term storage.
20X RT Enzyme Mix (50 × 20-µL reactions)	50 µL	500 µL	Store at -25°C to -15°C.

Method

1. Use up to 2 µg of total RNA per 20-µL reaction.
2. Allow the kit components to thaw on ice.
3. Referring to the table below, calculate the volume of components needed to prepare the required number of reactions.

Component	Volume per reaction	
	+RT reaction	- RT control
2X RT Buffer Mix	10.0 µL	10.0 µL
20X RT Enzyme Mix	1.0 µL	—
RNA sample	up to 9 µL	up to 9 µL
Nuclease-free H ₂ O	Q.S. ^[1] to 20 µL	Q.S. ^[1] to 20 µL
Total per reaction	20.0 µL	20.0 µL

^[1] Quantity Sufficient

4. Aliquot the RT reaction mix into a plate or tubes.
5. Seal the plates or tubes with the appropriate seal or caps.
6. Briefly centrifuge the plate or tubes to spin down the contents and to eliminate any air bubbles.
7. Place the plate or tubes on ice until you are ready to start the reverse transcription reaction.
8. Incubate the reaction for 37°C for 60 minutes. Stop the reaction by heating to 95°C for 5 minutes and hold at 4°C. For convenience, the incubation may be performed in a thermal cycler.
9. The cDNA is ready for use in real-time PCR applications or long-term storage in a freezer (-25°C to -15°C).

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.



Manufacturer: Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 | Vilnius, Lithuania

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.

Revision history: Revision history of Pub. No. 4387949

Revision	Date	Description
D	18 June 2018	Product line extension
C	21 March 2016	Format, style, and legal updates
B	15 August 2007	Baseline for this revision history

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

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.