TaqMan[®] Microbial Assays—single tube

Catalog Numbers A39420, A41332, A41333

Pub. No. MAN0018536 Rev. A.0

Note: For safety and biohazard guidelines, see the "Safety" appendix in the *TaqMan*[®] *Gene Expression Assays User Guide*—*single-tube assays* (Pub. No. 4333458). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

This document is intended as a benchtop reference for experienced users. See "Related documentation" on page 2 for resources that contain detailed instructions and troubleshooting.

Procedural guidelines

Guidelines for nucleic acid isolation

- For detailed instructions about nucleic acid isolation for respiratory tract microbiota experiments using the MagMAX[™] Viral/Pathogen Ultra Nucleic Acid Isolation Kit, see the appropriate resources that are listed in "Related documentation" on page 2.
- A 400-µL sample input yields approximately 50 µL of purified sample. Depending on the number of assays to be tested, multiple 400-µL sample aliquots may be needed.

Guidelines for 1-Step RT-PCR

- Use TaqPath[™] 1-Step RT-qPCR Master Mix, CG (Cat. No. A15299).
- Use purified, nondegraded total nucleic acid that is dissolved in PCR-compatible buffer.
- Ensure that the input nucleic acid is free of RNase activity and inhibitors of reverse transcription (RT) and PCR.
- To maximize sensitivity, prepare each reaction so that approximately half of the reaction volume is purified total nucleic acid.
- Protect the assays from light and store as indicated until ready for use. Excessive exposure to light can negatively affect the fluorescent probes of the assays.
- (*Optional*) Run technical replicates in triplicate to identify outliers.

Perform 1-Step RT-PCR

Prepare 1-Step RT-PCR reactions

Thaw the respiratory tract nucleic acid samples on ice. Resuspend the samples by inverting the tube, then gently vortexing.

- Gently mix the bottle of TaqPath[™] 1-Step RT-qPCR Master Mix, CG.
- 2. Combine the following components for the number of required reactions plus 10% overage:

IMPORTANT! For optimal results, prepare the reaction plate on ice.

	Volume for 1 reaction	
Component	Standard 96– well or 48– well Plates	384–well Plate or 96– well Fast Plate
TaqPath™ 1-Step RT-qPCR Master Mix, CG (4X)	5.0 µL	2.5 µL
TaqMan [®] Microbial Assay (20X)	1.0 µL	0.5 µL
Nuclease-free Water	4.0 µL	2.0 µL
Total PCR Reaction Mix volume	10.0 µL	5.0 µL

- 3. Vortex the PCR Reaction Mix, then centrifuge briefly.
- **4.** Combine the following components in each well of an optical reaction plate.

	Volume per well	
Component	Standard 96– well or 48– well Plates	384-well Plate or 96- well Fast Plate
PCR Reaction Mix (see step 2)	10.0 µL	5.0 µL
Extracted respiratory tract nucleic acid sample <i>or</i> elution buffer for NTC	10.0 µL	5.0 µL
Total volume	20.0 µL	10.0 µL

- 5. Seal the plate with a MicroAmp[™] Optical Adhesive Film, then vortex briefly to mix the contents.
- **6.** Centrifuge the plate briefly to collect the contents at the bottom of the wells.

IMPORTANT! Run the plate within 2 hours of preparation, or store the plate at 2–8°C for up to 24 hours.



Set up and run the real-time PCR instrument

See the appropriate instrument user guide for detailed instructions to program the thermal-cycling conditions or to run the plate.

Note: The instrument must be configured with the block appropriate for the plate type.

1. Select the Fast cycling mode.

IMPORTANT! The cycling mode depends on the master mix that is used in the reaction. The cycling mode does not depend on a Standard or a Fast plate format.

2. Set up the thermal protocol for your instrument.

Table 1 Thermal protocol for TaqPath[™] 1-Step RT-qPCR Master Mix, CG (StepOne[™], StepOnePlus[™], ViiA[™] 7, 7500 Real-Time PCR System, 7500 Fast Real-Time PCR System and QuantStudio[™] systems with fast cycling mode)

Stage	Step	Temperature	Time
Hold	UNG incubation ^[1]	25°C	2 minutes
Hold	Reverse transcription	50°C	15 minutes
Hold	Activation	95°C	2 minutes ^[2]
Cycling (40 cycles)	Denaturation	95°C	3 seconds
	Anneal/ Extension	60°C	1 minute

Heat-labile UNG is completely inactivated during the initial ramp to 95°C.
Required for RT inactivation, initial denaturation, and activation of the DNA polymerase.

- 3. Set the reaction volume appropriate for the reaction plate.
 - 96-well Standard (0.2-mL) Plate: 20 µL
 - 96-well Fast (0.1-mL) Plate and 384–well Plate: $10\ \mu L$
- 4. Load the plate into the real-time PCR instrument.
- 5. Start the run.

Analyze the results

For more information about analyzing the results, see the appropriate resources that are listed in "Related documentation" on page 2.

Use the absolute or relative quantification ($\Delta\Delta C_t$) methods (without target normalization) to analyze results.

The general guidelines for analysis include:

- View the amplification plot; then, if needed:
 - Adjust the baseline and threshold values.
 - Remove outliers from the analysis.
- In the well table or results table, view the C_t values for each well and for each replicate group, if applicable.

Related documentation

Document	Publication Number
Respiratory Tract Microbiota Profiling Experiments using OpenArray™ Application Guide	MAN0017952
Respiratory Tract Microbiota Profiling Experiments using TaqMan® Array Cards Application Guide	MAN0017951
TaqMan [®] Gene Expression Assays User Guide— single-tube assays	4333458
<i>Isolation of Nucleic Acid for Respiratory Tract</i> <i>Microbiota Profiling Experiments Quick Reference</i>	MAN00018526

Customer and technical support

Visit **thermofisher.com/support** for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation
 - User guides, manuals, and protocols



Life Technologies Corporation | 6055 Sunol Blvd | Pleasanton, CA 94566 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.

EARLY ACCESS DISCLAIMER: Customer acknowledges that this product is not a commercially released product and is still under development. THIS PRODUCT AND ASSOCIATED MATERIALS ARE SUPPLIED "AS IS" WITHOUT ANY WARRANTY, EXPRESS OR IMPLIED, AND ALL WARRANTIES ARE EXPRESSLY DISCLAIMED, INCLUDING WITHOUT LIMITATION THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, WHETHER ARISING FROM A STATUTE OR OTHERWISE IN LAW OR FROM A COURSE OF DEALING OR USAGE OF TRADE. Nothing herein shall be deemed to be or imply a duty for Life Technologies and/or its affiliate(s) to further develop this product or to release a commercial version thereof.

Revision history: Pub. No. MAN0018536

Revision	Date	Description
A.0	18 June 2019	New document.

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.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. TaqMan is a registered trademark of Roche Molecular Systems, Inc., used under permission and license.

- Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

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.

