


TrueMark™ Respiratory Panel 2.0 Amplification Control

Catalog Number A48101

Pub. No. MAN0019505 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

TrueMark™ Respiratory Panel 2.0 Amplification Control contains a linearized multi-target plasmid with target sequences for each available respiratory tract microbiota profiling assay, including SARS-CoV-2 virus assays. The plasmid also contains target sequences for TaqMan™ Universal RNA Spike In/Reverse Transcription (Xeno) Control, TaqMan™ Universal Extraction Control Organism (*B. atrophaeus*), and the human RNase P RPPH1 gene. The control can be included in respiratory tract microbiota profiling experiments as a positive control for panel-specific amplification.

TrueMark™ Respiratory Panel 2.0 Amplification Control is supplied at a concentration of 1×10^5 copies/μL. During real-time PCR, the control can be used as a stand-alone sample to verify assay performance and to help with troubleshooting.

See “Related documentation” on page 2 for resources that contain information about respiratory tract microbiota profiling workflows, including:

- A list of all available TaqMan™ assays for respiratory tract microbiota profiling
- Detailed instructions for respiratory tract microbiota profiling experiments using TaqMan™ Array Card, TaqMan™ OpenArray™ Plate, or single-tube assay formats
- Troubleshooting

Contents and storage

Component	Amount	Storage ^[1]
TrueMark™ Respiratory Panel 2.0 Amplification Control	250 μL (1×10^5 copies/μL)	-25°C to -15°C

^[1] Refer to label for expiration date.

Guidelines for use

- Before use, thaw the amplification control, vortex to thoroughly mix the contents, then centrifuge briefly to spin down the contents.
- Use the appropriate volume of the amplification control for your microbiota profiling experiment.

Format	Control volume
TaqMan™ Array Card	10.0 μL
Format 112 TaqMan™ OpenArray™ Plate	2.5 μL/well; 5.0 μL
Single-tube assay	96-well or 48-well standard plate: 14.0 μL 384-well plate or 96-well fast plate: 7.0 μL

- See “Related documentation” on page 2 for resources that include specific instructions for using the control.

Related documentation

Document	Pub. No.
<i>Respiratory Tract Microbiota Profiling Experiments v2 using TaqMan™ Array Cards Application Guide</i>	MAN0019507
<i>TaqMan™ Microbial Assays—single tube Quick Reference</i>	MAN0018536



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](https://www.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.

Revision history: Pub. No. MAN0019505

Revision	Date	Description
A.0	29 October 2020	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.

©2020 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.