

# TrueMark™ Comprehensive Microbiota Control

Catalog Numbers A50382 and A50383

Pub. No. MAN0024964 Rev. E



**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).

## Product description

The Applied Biosystems™ TrueMark™ Comprehensive Microbiota Control contains a linearized multi-target plasmid pool with microbial target sequences commonly used on TrueMark™ OpenArray™ panels and TaqMan™ Array Card panels.

The control has targets relevant to many microbial profiling applications, including respiratory tract, urinary tract, vaginal microbiota, and antibiotic resistance research.

For a list of targets, go to: <https://www.thermofisher.com/order/catalog/product/A50382>.

The control can be included as a positive control for panel-specific amplification. The high-concentration control (Cat. No. A50383) can also be used to determine parameters necessary for analytical confirmation (e.g. limit of detection and  $C_t$  cut-off values).

The plasmid also contains target sequences for the TrueMark™ Universal RNA Spike In/Reverse Transcription (Xeno) Control, TrueMark™ Universal Extraction Control Organism (*B. atrophaeus*), and the human RNase P RPPH1 gene. During real-time PCR, the control can be used as a stand-alone sample to verify analytical performance of an assay and to help with troubleshooting.

This product is updated on a regular basis to support coverage of the most popular microbial targets. Analytical performance is measured to support consistency between versions. Version 1 of this product was first available in April 2021 and contained 471 microbial assay targets. Version 2 became available in October 2021 and contained an additional 83 microbial assay targets. Version 3 became available in March 2023 and contained an additional 139 microbial and antibiotic resistant targets. Version 4 became available in February 2025 and contains an additional 150 microbial targets.

## Contents and storage

Store the TrueMark™ Comprehensive Microbiota Control at –25°C to –15°C.

Name	Cat. No.	Concentration	Volume
TrueMark™ Comprehensive Microbiota Control	A50382	$1 \times 10^5$ copies/μL	1,000 μL
	A50383	$5 \times 10^7$ copies/μL	50 μL

## Guidelines

- Products are stable for two years from date of manufacture when stored at –20°C.
- Before first use, thaw the control, vortex to thoroughly mix the contents, then centrifuge briefly to collect the contents at the bottom of the tube
- Create aliquots to reduce the number of freeze-thaw cycles.
- Do not exceed three freeze-thaw cycles.
- Aliquots are stable for four months when stored at 4°C, if the products are still within two years of the manufacture date.
- Use the appropriate volume of the control at  $1 \times 10^5$  copies/μL for the microbiota profiling experiment.

Format	Volume of control
TaqMan™ Array Card	10.0 μL
TrueMark™ OpenArray™ Plate	1.0 μL

- For more information about using the control in your microbiota profiling experiment, see “Related documentation” on page 2.

## Related documentation

Document	Publication number
<i>Respiratory Tract Microbiota Profiling Experiments using TaqMan™ Array Cards Application Guide</i>	MAN0017951
<i>Respiratory Tract Microbiota Profiling Experiments v2 using OpenArray™ Plates Application Guide</i>	MAN0019506
<i>Respiratory Tract Microbiota Profiling Experiments v2 using TaqMan™ Array Cards Application Guide</i>	MAN0019507
<i>TrueMark™ Controls with TaqMan™ Array Cards Quick Reference</i>	MAN0025035
<i>TrueMark™ Controls with TrueMark™ OpenArray™ Plates Quick Reference</i>	MAN0025036
<i>TrueMark™ OpenArray™ Plates for Respiratory Tract Microbiota Profiling Experiments User Guide</i>	MAN0017952
<i>TrueMark™ Pre-Designed and Custom Multiplex Panels User Guide</i>	MAN1000315
<i>Urinary Tract Microbiota Profiling Experiments Application Guide</i>	MAN0017750
<i>Vaginal Microbiota Profiling Experiments Application Guide</i>	MAN0015669

## Limited product warranty

Life Technologies Corporation and its affiliates 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](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have questions, contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



Thermo Fisher Scientific GENEART GmbH | Im Gewerbepark B35 | 93059 Regensburg | Germany

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](http://thermofisher.com/symbols-definition).

Revision history: Pub. No. MAN0024964 E

Revision	Date	Description
E	16 April 2025	<ul style="list-style-type: none"><li>Information for Version 3 and Version 4 updates was added to "Product description".</li><li>"Related documentation" was updated.</li></ul>
D.0	27 October 2023	<ul style="list-style-type: none"><li>TaqMan™ branding was updated to TrueMark™ throughout document.</li><li>Shelf life was updated to two years in "Guidelines" on page 1.</li></ul>
C.0	24 September 2021	Updated the product description to include an annual review of the microbial targets.
B.0	6 May 2021	Quick reference documents added.
A.0	12 March 2021	New document for the TrueMark™ Comprehensive Microbiota Control.

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 these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2021-2025 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 trademark of Roche Molecular Systems, Inc., used under permission and license.