


# TaqMan® Urinary Tract Microbiota Amplification Control

Catalog Number A39174

Pub. No. MAN0017753 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](http://thermofisher.com/support).

## Product information

The TaqMan® Urinary Tract Microbiota Amplification Control (Cat. No. A39174) contains a linearized multi-target plasmid with target sequences for each available urinary tract microbiota profiling assay. See the *Urinary Tract Microbiota Profiling Experiments Application Guide* (Pub. No. MAN0017750) for the full list of available TaqMan® urinary tract profiling assays. The plasmid also contains target sequences for Xeno DNA and human RNase P RPPH1 genes, for a general control for the sample preparation process.

The TaqMan® Urinary Tract Microbiota Amplification Control can be included in urinary tract microbiota profiling experiments to verify assay performance and to help with troubleshooting.

The TaqMan® Urinary Tract Microbiota Amplification Control is supplied at a plasmid concentration of 10<sup>5</sup> copies/μL. The amplification control is designed for use with urinary tract microbiota profiling protocols, see “Related documentation” on page 2.

See “Related documentation” on page 2 for resources that contain detailed instructions and troubleshooting for urinary tract microbiota profiling OpenArray™ experiments.

## Contents and storage

Component	Amount	Storage
TaqMan® Urinary Tract Microbiota Amplification Control	5 × 10 μL (sufficient for up to 20 reactions)	-30°C to -15°C (long-term) 4°C (up to 2 months)

## Guidelines for use

These guidelines apply to all OpenArray™ plate formats available for urinary tract microbiota profiling experiments.

- Before use, thaw the amplification control, vortex to thoroughly mix the contents, then centrifuge briefly to spin down the contents.
- Use the amplification control at the concentration supplied; no dilution is required.
- Use TaqMan® OpenArray™ Real-Time PCR Master Mix for urinary tract microbiota profiling experiments.
- Add 2.5 μL of amplification control and 2.5 μL of master mix to each Amplification Control well of the OpenArray™ 384-well Sample Plate as designated in your sample layout.
- Add the DNA samples and master mix to the appropriate wells of the same OpenArray™ 384-well Sample Plate, then continue the OpenArray™ experiment protocol as described in the *Urinary Tract Microbiota Profiling Experiments Application Guide* (Pub. No. MAN0017750).

## Related documentation

Document	Pub. No.
<i>Urinary Tract Microbiota Profiling Experiments Application Guide</i>	MAN0017750
<i>DNA isolation for Urinary Tract Microbiota Profiling Experiments Quick Reference</i>	MAN0017751
<i>OpenArray™ Urinary Tract Microbiota Profiling Experiments Quick Reference</i>	MAN0017752
<i>QuantStudio™ 12K Flex Real-Time PCR System: OpenArray™ Experiments User Guide</i>	4470935
<i>OpenArray™ Sample Tracker Software Quick Reference</i>	4460657
<i>OpenArray™ AccuFill™ System User Guide</i>	4456986



**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:** Pub. No. MAN0017753

Revision	Date	Description
A.0	06 August 2018	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.

©2018 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.