


# TaqMan™ Custom DNA Controls

Pub. No. MAN0024963 Rev. B.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 description

The Applied Biosystems™ TaqMan™ Custom DNA Controls contain a linearized multi-target plasmid pool used for microbial detection with TaqMan™ Array Cards or TaqMan™ OpenArray™ Plates.

The controls can be included as a positive control for panel-specific amplification. The high concentration ( $5 \times 10^7$  copies/μL) products can also be used to determine the parameters necessary for analytical confirmation (e.g. limit of detection and  $C_t$  cut-off values).

During real-time PCR, the control can be used as a stand-alone sample to verify assay performance and to help with troubleshooting.

## Contents and storage

Store the TaqMan™ Custom DNA Controls at  $-25^\circ\text{C}$  to  $-15^\circ\text{C}$ .

| Name                       | Cat. No. | Number of targets | Concentration             | Volume   |
|----------------------------|----------|-------------------|---------------------------|----------|
| TaqMan™ Custom DNA Control | A50319   | $\leq 20$         | $1 \times 10^5$ copies/μL | 1,000 μL |
|                            | A50375   | 21–30             | $1 \times 10^5$ copies/μL | 1,000 μL |
|                            | A50376   | 31–50             | $1 \times 10^5$ copies/μL | 1,000 μL |
|                            | A50377   | $> 50$            | $1 \times 10^5$ copies/μL | 1,000 μL |
|                            | A50378   | $\leq 20$         | $5 \times 10^7$ copies/μL | 50 μL    |
|                            | A50379   | 21–30             | $5 \times 10^7$ copies/μL | 50 μL    |
|                            | A50380   | 31–50             | $5 \times 10^7$ copies/μL | 50 μL    |
|                            | A50381   | $> 50$            | $5 \times 10^7$ copies/μL | 50 μL    |

## Guidelines

- Products are stable for one year from date of manufacture when stored at  $-20^\circ\text{C}$ .
- 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  $4^\circ\text{C}$ , if the products are still within one year 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           |
| TaqMan™ 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>TaqMan™ Controls with TaqMan™ Array Cards Quick Reference</i>   | MAN0025035         |
| <i>TaqMan™ Controls with TaqMan™ OpenArray™ Plates Quick Reference</i>                                   | MAN0025036         |
| <i>Respiratory Tract Microbiota Profiling Experiments using TaqMan™ Array Cards Application Guide</i>    | MAN0017951         |
| <i>Respiratory Tract Microbiota Profiling Experiments v2 using TaqMan™ Array Cards Application Guide</i> | MAN0019507         |
| <i>Respiratory Tract Microbiota Profiling Experiments using OpenArray™ Application Guide</i>             | MAN0017952         |
| <i>Respiratory Tract Microbiota Profiling Experiments v2 using OpenArray™ Application Guide</i>          | MAN0019506         |
| <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/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](http://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](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).

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. MAN0024963

| Revision | Date          | Description                                       |
|----------|---------------|---|
| B.0      | 6 May 2021    | Quick reference documents added.                  |
| A.0      | 12 March 2021 | New document for the TaqMan™ Custom DNA Controls. |

**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 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.