


# TaqMan™ Universal DNA Spike In Control

Universal control for DNA extraction and qPCR workflows

Catalog Number A39175

Pub. No. MAN0017852 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](https://www.thermofisher.com/support).

## Contents and storage

Contents	Amount	Storage
TaqMan™ Universal DNA Spike In Control	1 mL 2 × 10 <sup>5</sup> copies/μL	-25°C to -15°C

## Product information

The TaqMan™ Universal DNA Spike In Control serves as an exogenous process control of known concentration that can be added to a sample at various stages throughout the qPCR workflow. Coupled with the proprietary TaqMan™ Assay for the Xeno™ DNA control (Assay ID Ac00010014\_a1), the TaqMan™ Universal DNA Spike In Control can be used to monitor the recovery efficiency for the DNA extraction and purification process. The control can also be used to indicate the presence of PCR inhibitors in molecular detection workflows. This level of verification is of particular importance when working with sample types that tend to have higher levels of PCR inhibitors. This verification layer helps to ensure that the PCR results are accurate, and it reduces the likelihood of false negatives.

The TaqMan™ Universal DNA Spike In Control sequence does not overlap with any known microbial organism sequence, enabling its use with any microbial PCR workflow. This product contains a sequence for AmpC beta-lactamase and may amplify for any assays designed to detect this antibiotic resistance target. For example, TaqMan™ Assay ID Ba04646128\_s1 will amplify in any sample that contains the TaqMan™ Universal DNA Spike In Control. Typically, the control is added during DNA isolation, then carried through the remainder of the qPCR workflow. The ready-to-use control is compatible with any target assay, master mix, or sample preparation reagents.

## Guidelines for use

- TaqMan™ Universal DNA Spike In Control is supplied ready to use.
- Add 10 μL of the control (2 × 10<sup>6</sup> copies) per reaction to the lysis solution used for DNA isolation. Continue the DNA isolation procedure without further changes.
- Use in conjunction with the TaqMan™ Assay for the Xeno™ DNA control (Assay ID Ac00010014\_a1).
- The TaqMan™ Universal DNA Spike In Control has an expected C<sub>t</sub> range of 19-23.

## 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](https://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](https://www.thermofisher.com/support).

[thermofisher.com/support](https://www.thermofisher.com/support) | [thermofisher.com/askaquestion](https://www.thermofisher.com/askaquestion)



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. MAN0017852

Revision	Date	Description
B.0	29 March 2021	Updated product information.
A.0	6 August 2018	New document.

**Important Licensing Information:** This product may be covered by one or more Limited Use Label Licenses. By use of this product, 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.

**For Research Use Only. Not for use in diagnostic procedures.**