

TaqMan® Fast Universal PCR Master Mix, no AmpErase™ UNG

Catalog Numbers 4352042, 4364103, 4366072, 4366073, and 4367846

Pub. No. 4357649 Rev. E

Note: For safety and biohazard guidelines, see the “Safety” appendix in the *TaqMan® Fast Universal PCR Master Mix, no AmpErase™ UNG User Guide* (Pub. No. 4351891). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Contents and storage

Cat. No.	Number of reactions	Contents	Storage
4352042	250	2 × 1.25 mL	2–8°C ^[1]
4364103	2,500	20 × 1.25 mL	
4366072	500	4 × 1.25 mL	
4366073	1,250	10 × 1.25 mL	
4367846	5,000	50 mL	

^[1] See packaging for expiration date.

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 found on Life Technologies' website at 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.



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. 4357649

Revision	Date	Description
E	04 September 2018	Updated manufacturing address, branding, licensing, trademarks, general style and format.
D	July 2011	Baseline for this revision.

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.