

TaqMan® Universal Master Mix II

Catalog Numbers 4440040, 4440041, 4440043, 4440047, 4440048, 4440049, 4440038, 4440039, 4440042, 4440044, 4440045, 4440046

Pub. No. 4428175 **Rev.** C

Note: For safety and biohazard guidelines, see the “Safety” appendix in the *TaqMan® Universal Master Mix II User Guide* (Pub. No. 4428173). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Product description

Applied Biosystems™ TaqMan® Universal Master Mix II is a mix of components necessary to perform real-time polymerase chain reaction (PCR). TaqMan® Universal Master Mix II can be used to amplify complementary DNA (cDNA) and DNA targets for quantitation, genotyping, and other applications. The mix is available with or without UNG. TaqMan® Universal Master Mix II is optimized for use with TaqMan® primers and TaqMan® probes designed according to Applied Biosystems™ guidelines. It can be used with custom TaqMan® assays and predesigned assays including TaqMan® Gene Expression Assays, TaqMan® MicroRNA Assays, TaqMan® Drug Metabolism Genotyping Assays and TaqMan® SNP Genotyping Assays.

Contents and storage

Item	Cat. No.	Amount	Storage ^[1]
TaqMan® Universal Master Mix II, with UNG ^[2]	4440042	1 × 1 mL	2–8°C
	4440038	1 × 5 mL	
	4440044	2 × 5 mL	
	4440045	5 × 5 mL	
	4440046	10 × 5 mL	
	4440039	1 × 50 mL	
TaqMan® Universal Master Mix II, no UNG	4440043	1 × 1 mL	
	4440040	1 × 5 mL	
	4440047	2 × 5 mL	
	4440048	5 × 5 mL	
	4440049	10 × 5 mL	
	4440041	1 × 50 mL	

^[1] See packaging for expiration date.

^[2] Uracil-N-Glycosylase

Additional information

For detailed procedures, safety guidelines and component descriptions, see *TaqMan® Universal Master Mix II User Guide* (Pub. No. 4428173).

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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



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. 4428175

Revision	Date	Description
C	15 October 2018	Updated for manufacturer, general style, formatting, and branding.
B	July 2009	Baseline for this revision history.

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.