

TaqMan® OpenArray™ Genotyping Master Mix

Catalog Numbers 4404846

Pub. No. 4401098 Rev. E

Note: For safety and biohazard guidelines, see the “Safety” appendix in the *TaqMan® OpenArray™ Genotyping Getting Started Guide* (Pub. no. 4377476). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Contents and storage

Contents	Amount	Storage
TaqMan® OpenArray™ Genotyping Master Mix, 2X	5 mL	2–8°C

Product description

The 2X TaqMan® OpenArray™ Genotyping Master Mix is optimized for SNP genotyping using TaqMan® OpenArray™ Real-Time PCR Plate with Genotyping Assays (Cat. No. 4471113).

Recommended genotyping PCR procedures

For detailed PCR reaction mix setup and PCR amplification parameters, see *TaqMan® OpenArray™ Genotyping Getting Started Guide* (Pub. No. 4377476).

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. 4401098

Revision	Date	Description
E	23 August 2018	Changed manufacturer address to Vilnius.
D	10 October 2017	Updated and rebranded.
C	December 2010	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.