

TrueMark™ Respiratory Panel 2.0, OpenArray™ Plate

Catalog Number A49044

Pub. No. MAN0019510 Rev. A.0

Note: For safety and biohazard guidelines, see the “Safety” appendix in the *Respiratory Tract Microbiota Profiling Experiments v2 using OpenArray™ Application Guide* (Pub. No. MAN0019506). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

For resources that contain detailed information about using the TrueMark™ Respiratory Panel 2.0, OpenArray™ Plate, see “Related documentation” on page 2.

Product description

TrueMark™ Respiratory Panel 2.0, OpenArray™ Plate (Cat. No. [A49044](#)) is an efficient, easy-to-use fixed-content OpenArray™ plate for the characterization of key respiratory tract microbial targets. The plate includes 35 TaqMan™ assays that have been optimized for detection of 32 respiratory tract viral and bacterial nucleic acid, including two assays that target SARS-CoV-2. The plate also includes control assays for TaqMan™ Universal Extraction Control Organism (*B. atrophaeus*), TaqMan™ Universal RNA Spike In/Reverse Transcription (Xeno) Control, and the human RNase P RPPH1 gene. For a complete list of assays included in the plate, see “TaqMan™ assays included in the TrueMark™ Respiratory Panel 2.0, OpenArray™ Plate” on page 1.

The assays perform well with total nucleic acid that is isolated from nasopharyngeal swab, nasopharyngeal aspirate, and bronchoalveolar lavage (BAL) research samples using the MagMAX™ Viral/Pathogen Ultra Nucleic Acid Isolation Kit.

TaqMan™ assay designs and assay target sequences have undergone rigorous bioinformatics selection and analysis to maximize strain coverage and minimize potential for off-target cross-reactivity. Qualified TaqMan™ assays have undergone performance testing to ensure that results are accurate with high levels of sensitivity and specificity.

Contents and storage

Table 1 TrueMark™ Respiratory Panel 2.0, OpenArray™ Plate (Cat. No. [A49044](#); format 112 assays and up to 24 samples)

Component	Amount	Storage
TrueMark™ Respiratory Panel 2.0, OpenArray™ Plate	1 plate	-25°C to -15°C

TaqMan™ assays included in the TrueMark™ Respiratory Panel 2.0, OpenArray™ Plate

The following assays are included in the TrueMark™ Respiratory Panel 2.0, OpenArray™ Plate.

Note: This format 112 TaqMan™ OpenArray™ Plate contains 35 TaqMan™ assays for 32 microbial targets, and the TaqMan™ Universal RNA Spike In/Reverse Transcription (Xeno) Control, plated in triplicate, and assays for the TaqMan™ Universal Extraction Control Organism (*B. atrophaeus*) and human RNase P RPPH1 gene controls plated in duplicate. Each TaqMan™ OpenArray™ Plate can be used for running up to 23 samples and one control sample.

Table 2 Assays for respiratory tract microbiota

Target organism	Assay name	Nucleic acid type	Assay ID
Bacteria			
<i>Bordetella bronchiseptica</i> / <i>parapertussis</i> / <i>pertussis</i>	Bordetella	DNA	Ba06439624_s1
<i>Bordetella pertussis</i>	B.pertussis	DNA	Ba06439623_s1
<i>Chlamydomphila pneumoniae</i>	C.pneumoniae	DNA	Ba06439616_s1
<i>Haemophilus influenzae</i>	H.influenzae	DNA	Ba06439625_s1
<i>Klebsiella pneumoniae</i> complex	K. pneumoniae	DNA	Ba04932083_s1
<i>Legionella pneumophila</i>	L.pneumophila	DNA	Ba06439617_s1
<i>Mycoplasma pneumoniae</i>	M.pneumoniae	DNA	Ba06439620_s1
<i>Staphylococcus aureus</i>	S.aureus	DNA	Ba04646259_s1
<i>Streptococcus pneumoniae</i>	S.pneumoniae	DNA	Ba06439619_s1

Target organism	Assay name	Nucleic acid type	Assay ID
Virus			
Adenovirus	AdV_1of2	DNA	Vi99990001_po
Adenovirus	AdV_2of2	DNA	Vi99990002_po
Human Coronavirus 229E	CoV_229E	RNA	Vi06439671_s1
Human Coronavirus HKU1	CoV_HKU1	RNA	Vi06439674_s1
Human Coronavirus NL63	CoV_NL63	RNA	Vi06439673_s1
Human Coronavirus OC43	CoV_OC43	RNA	Vi06439646_s1
Human Enterovirus (pan assay)	EV_pan	RNA	Vi06439631_s1
Human Enterovirus D68	EV_D68	RNA	Vi06439669_s1
Human Metapneumovirus (hMPV)	hMPV	RNA	Vi99990004_po
Human Parainfluenza virus 1	hPIV1	RNA	Vi06439642_s1
Human Parainfluenza virus 2	hPIV2	RNA	Vi06439672_s1
Human Parainfluenza virus 3	hPIV3	RNA	Vi06439670_s1
Human Parainfluenza virus 4	hPIV4	RNA	Vi99990005_po
Human Respiratory Syncytial Virus A (RSVA)	RSVA	RNA	Vi99990014_po
Human Respiratory Syncytial Virus B (RSVB)	RSVB	RNA	Vi99990015_po
Human Rhinovirus 1/2	RV_1of2	RNA	Vi99990016_po
Human Rhinovirus 2/2	RV_2of2	RNA	Vi99990017_po
Human herpesvirus 4 (HHV4 – Epstein-Barr Virus)	HHV4	DNA	Vi06439675_s1
Human herpesvirus 5 (HHV5 – Cytomegalovirus)	HHV5	DNA	Vi06439643_s1
Human herpesvirus 6 (HHV6)	HHV6	DNA	Vi06439627_s1
Influenza A	Flu_A_pan	RNA	Vi99990011_po
Influenza A/H1-2009	Flu_A_H1	RNA	Vi99990009_po
Influenza A/H3	Flu_A_H3	RNA	Vi99990010_po
Influenza B	Flu_B_pan	RNA	Vi99990012_po
Severe Acute Respiratory Syndrome coronavirus 2 (SARS-CoV-2); S protein gene	SARS-CoV2S	RNA	Vi07918636_s1
Severe Acute Respiratory Syndrome coronavirus 2 (SARS-CoV-2); N protein gene	SARS-CoV2N	RNA	Vi07918637_s1

Table 3 Assays for respiratory tract microbiota controls

Control name	Assay name	Nucleic acid type	Assay ID
TaqMan™ Universal Extraction Control Organism (<i>B. atrophaeus</i>)	B.atrophaeus	DNA	Ba06596576_s1
TaqMan™ Universal RNA Spike In/Reverse Transcription (Xeno) Control	Xeno	RNA	Ac00010014_a1
Human RNase P RPPH1 gene	RPPH1	DNA	Hs04930436_g1

Related documentation

Document	Publication Number
<i>Respiratory Tract Microbiota Profiling Experiments v2 using OpenArray™ Application Guide</i>	MAN0019506
<i>Respiratory Tract Microbiota Profiling Experiments v2 using OpenArray™ Quick Reference</i>	MAN0019509
<i>Isolation of Nucleic Acid for Respiratory Tract Microbiota Profiling Experiments Quick Reference</i>	MAN0018526

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.



Life Technologies Corporation | 6055 Sunol Blvd | Pleasanton, CA 94566 USA

For descriptions of symbols on product labels or product documents, go to 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. MAN0019510

Revision	Date	Description
A.0	5 January 2021	New document.

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.