

# TaqMan® Array Respiratory Tract Microbiota Comprehensive Card

Catalog Number A41238

Pub. No. MAN0018632 Rev. A.0

**Note:** For safety and biohazard guidelines, see the “Safety” appendix in the *Respiratory Tract Microbiota Profiling Experiments using TaqMan® Array Cards Application Guide* (Pub. No. MAN0017951). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

## Product description

TaqMan® Array Respiratory Tract Microbiota Comprehensive Card is an efficient, easy-to-use TaqMan® Array Card for the characterization of key respiratory tract microbes. The array card includes TaqMan® assays that have been optimized for detection of 42 respiratory tract viral, bacterial, and fungal microbes. The array card also includes control assays for TaqMan® Universal Extraction Control Organism (*B. atrophaeus*), TaqMan® Universal RNA Spike In/Reverse Transcription (Xeno) Control, the human RNase P RPPH1 gene, and the human 18S ribosomal RNA gene. For a complete list of assays included in the TaqMan® Array Respiratory Tract Microbiota Comprehensive Card, see Table 1.

The assays perform well with total nucleic acid that is isolated from nasopharyngeal swab, nasopharyngeal aspirate, and bronchoalveolar lavage (BAL) samples using the MagMAX™ Viral/Pathogen Ultra Nucleic Acid Isolation Kit (Cat. No. A42356).

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 also undergo performance testing to ensure that results are accurate and reproducible with high levels of sensitivity and specificity.

See “Related documentation” on page 2 for resources that contain detailed information about using the TaqMan® Array Respiratory Tract Microbiota Comprehensive Card.

## Contents and storage

Component	Amount	Storage
TaqMan® Array Respiratory Tract Microbiota Comprehensive Card	10 cards <sup>[1]</sup>	2°C to 8°C

<sup>[1]</sup> Minimum order.

## TaqMan® assays for respiratory tract microbiota profiling

The following assays are included in the TaqMan® Array Respiratory Tract Microbiota Comprehensive Card.

**Table 1** Assays for respiratory tract microbiota targets

Target organism	Assay name	Nucleic acid type	Assay ID
<b>Bacteria</b>			
<i>Bordetella bronchiseptica</i> / <i>parapertussis</i> / <i>pertussis</i>	Bordetella	DNA	Ba06439624_s1
<i>Bordetella holmesii</i>	B.holmesii	DNA	Ba06439621_s1
<i>Bordetella pertussis</i>	B.pertussis	DNA	Ba06439623_s1
<i>Chlamydophila pneumoniae</i>	C.pneumoniae	DNA	Ba06439616_s1
<i>Coxiella burnetii</i>	C.burnetii	DNA	Ba06439618_s1
<i>Haemophilus influenzae</i>	H.influenzae	DNA	Ba06439625_s1
<i>Klebsiella pneumoniae</i>	K.pneumoniae	DNA	Ba04932083_s1
<i>Legionella pneumophila</i>	L.pneumophila	DNA	Ba06439617_s1
<i>Moraxella catarrhalis</i>	M.catarrhalis	DNA	Ba06439622_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
<b>Fungus</b>			
<i>Pneumocystis jirovecii</i>	P.jirovecii	DNA	Fn06439626_s1

Target organism	Assay name	Nucleic acid type	Assay ID
<b>Virus</b>			
Adenovirus	AdV_1of2	DNA	Vi99990001_po
Adenovirus	AdV_2of2	DNA	Vi99990002_po
Human Bocavirus	HBoV	DNA	Vi99990003_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 Parechovirus	HPeV	RNA	Vi99990006_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 3 (HHV3 – Varicella zoster Virus)	HHV3	DNA	Vi06439647_s1
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
Measles virus	Measles	RNA	Vi99990013_po
Middle East Respiratory Syndrome coronavirus (MERS)	MERS_CoV	RNA	Vi06439644_s1
Mumps virus	Mumps	RNA	Vi06439657_s1
Severe Acute Respiratory Syndrome coronavirus (SARS)	SARS_CoV	RNA	Vi06439634_s1

**Table 2** 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
Human 18S ribosomal RNA gene <sup>[1]</sup>	18S	DNA	Hs99999901_s1

<sup>[1]</sup> Human 18S ribosomal RNA is used as a manufacturing control for the array card and is not an assay specific to Respiratory Tract Microbiota profiling.

## Related documentation

Document	Pub. No.
<i>Respiratory Tract Microbiota Profiling Experiments using TaqMan® Array Cards Application Guide</i>	MAN0017951
<i>Respiratory Tract Microbiota Profiling Experiments using TaqMan® Array Cards Quick Reference</i>	MAN0018528
<i>Isolation of Nucleic Acid for Respiratory Tract Microbiota Profiling Experiments Quick Reference</i>	MAN00018526

## 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](http://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](http://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. MAN0018632

Revision	Date	Description
A.0	27 August 2019	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.

©2019 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.