


TaqMan™ 2019-nCoV Control Kit v1

For use with TaqMan™ 2019-nCoV Assay Kit v1

Catalog Number A47533

Pub. No. MAN0019097 Rev. A.0

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product description

Applied Biosystems™ TaqMan™ 2019-nCoV Control Kit v1 (Cat. No. A47533) is a synthetic positive control that contains target sequences for each of the assays included in the TaqMan™ 2019-nCoV Assay Kit v1 (Cat. No. A47532).

- The control includes synthetic DNA target sequences for three SARS-CoV-2 genes (ORF1ab, S protein, and N protein) and the Human RNase P RPPH1 gene.
- During RT-PCR, use 1 µL of 2019-nCoV Control v1 as the sample for each positive-control reaction.
- The control can be used if needed to verify assay performance and to help with troubleshooting.

For more information about genetic analysis solutions available for SARS-CoV-2, go to thermofisher.com/coronavirus.

Contents and storage

Contents	Target Sequence	Amount ^[1]	Concentration	Storage
2019-nCoV Control v1	ORF1ab	50 µL	1 × 10 ⁴ copies/µL	-30°C to -10°C
	S protein			
	N protein			
	Human RNase P RPPH1			

^[1] Sufficient for 50 positive-control reactions. We recommend 3 positive control reactions per plate, one for each 2019-nCoV assay.

Guidelines for use

- Include the RNase P assay as an internal positive control for each of the positive-control reactions.
- We recommend that you include three positive-control reactions per plate, one for each of the 2019-nCoV assays.

Table 1 Recommended positive-control reactions (per plate)

Positive-control reaction	Volume of 2019-nCoV Control v1 per reaction
1. 2019-nCoV (ORF1ab) assay + RNase P assay	1 µL
2. 2019-nCoV (S Protein) assay + RNase P assay	1 µL
3. 2019-nCoV (N Protein) assay + RNase P assay	1 µL
Total volume of 2019-nCoV Control v1 required per plate	3 µL

- For guidelines about interpreting the results of the positive-control reactions, see *TaqMan™ 2019-nCoV Assay Kit v1 Product Information Sheet* (Pub. No. MAN0019096).

Related documentation

Document	Publication Number
<i>TaqMan™ 2019-nCoV Assay Kit v1 Product Information Sheet</i>	MAN0019096
<i>TaqMan™ Gene Expression Assays User Guide—single-tube assays</i>	4333458

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

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. MAN0019097

Revision	Date	Description
A.0	2 March 2020	New document.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.