

# Real-Time PCR Detection of SARS-CoV-2 on Food Packaging and Environmental Surfaces

## USER GUIDE

for use with:

PrepSEQ™ Nucleic Acid Extraction Kit for Food and Environmental Testing

KingFisher™ Flex-96 Deep-Well Magnetic Particle Processor

MagMAX™ Express-96 Deep Well Magnetic Particle Processor

TaqMan™ 2019-nCoV Assay Kit v1

TaqMan™ 2019-nCoV Control Kit v1

RNA UltraSense™ One-Step qPCR Master Mix

Applied Biosystems™ QuantStudio™ Design and Analysis Software v1.5.1 or later

Applied Biosystems™ 7500 Software SDS v.1.4.2 or later

Publication Number MAN0019499

Revision C.0



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](https://www.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. MAN0019499

Revision	Date	Description
C.0	12 October 2020	Addition of RNA UltraSense™ One-Step qPCR Master Mix
B.0	02 October 2020	Correction of sample volume on page 15
A.0	11 September 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.

**Trademarks:** 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. Thermomixer is a trademark of Eppendorf.

©2020 Thermo Fisher Scientific Inc. All rights reserved.

# Contents

■	<b>CHAPTER 1</b>	<b>Product information</b> .....	4
		Product description .....	4
		Required materials .....	5
		Materials for sample preparation and nucleic acid extraction .....	5
		Materials for PCR detection and analysis .....	6
		Procedural guidelines .....	8
		Guidelines for sample collection and storage .....	8
		Guidelines for the 2019-nCoV assay .....	8
		Workflow overview .....	9
■	<b>CHAPTER 2</b>	<b>Methods</b> .....	10
		Prepare the samples and extract RNA using PrepSEQ™ Nucleic Acid Extraction Kit for Food and Environmental Testing .....	10
		Perform 1-Step RT-PCR .....	12
		Analyze data .....	13
		Review results .....	14
		Troubleshooting .....	16
■	<b>APPENDIX A</b>	<b>Supplemental information</b> .....	18
		Good laboratory practices for PCR .....	18
		Symbol definitions .....	19
■	<b>APPENDIX B</b>	<b>Safety</b> .....	20
		Chemical safety .....	21
		Biological hazard safety .....	22
■	<b>APPENDIX C</b>	<b>Documentation and support</b> .....	23
		Food Safety support .....	23
		Related documentation .....	23
		Reference .....	23



# Product information

---

**IMPORTANT!** Before using this product, read and understand the information in the “Safety” appendix in this document.

---

## Product description

This guide describes the following *Performance Tested* workflow for detection of SARS-CoV-2 in food samples and environmental surfaces with the Applied Biosystems™ TaqMan™ 2019-nCoV Assay Kit v1 (Cat. No. A47532). This kit includes three TaqMan™ RT-PCR assays that target SARS-CoV-2 genes, and one positive control assay that targets the Human RNase P RPPH1 gene:

- Assays target three different viral genomic regions, reducing the risk of false negatives.
- Assays have undergone bioinformatic selection and analysis to specifically target sequences that are unique to SARS-CoV-2.
- The RNase P assay is run in duplex with the combined 2019-nCoV assays as an internal positive control.

The kit is used with the TaqMan™ 2019-nCoV Control Kit v1 (Cat. No. A47533) to monitor assay-specific amplification.

For more information about genetic analysis solutions available for SARS-CoV-2, go to [thermofisher.com/coronavirus](https://thermofisher.com/coronavirus).

## Required materials

Unless otherwise indicated, all materials are available through the Thermo Fisher Microbiology ordering process or [thermofisher.com](https://www.thermofisher.com). MLS: Fisher Scientific ([fisherscientific.com](https://www.fisherscientific.com)) or other major laboratory supplier.

**Note:** Parts may ship separately depending on configuration and storage conditions.

## Materials for sample preparation and nucleic acid extraction

**Table 1** PrepSEQ™ Nucleic Acid Extraction Kit

Contents	Cat. No. 4480466 (100 reactions)	Cat. No. 4428176 (300 reactions)	Storage <sup>[1]</sup>
Lysis Buffer	2 × 50 mL	6 × 50 mL	15°C to 30°C
Magnetic Particles	2 × 1.5 mL	6 × 1.5 mL	
Binding Solution (Isopropanol) <sup>[2]</sup>	1 empty bottle	3 empty bottles	
Wash Buffer Concentrate <sup>[3]</sup>	2 × 26 mL	6 × 26 mL	
Elution Buffer	25 mL	3 × 25 mL	
Proteinase K (PK) Buffer	50 mL	3 × 50 mL	-25°C to -15°C
Proteinase K, 20 mg/mL	1.25 mL	3 × 1.25 mL	

<sup>[1]</sup> Refer to the product label for the expiration date.

<sup>[2]</sup> Add ~35 mL of 100% isopropanol to the empty bottle before use.

<sup>[3]</sup> Add 74 mL of 95% ethanol before use.

**Table 2** Magnetic particle processor

Item	Source
<b>KingFisher™ Flex-96 instrument and accessories</b>	
KingFisher™ Flex-96 Deep-Well Magnetic Particle Processor	A32681
KingFisher™ Flex -96 Deep-Well Heating Block	<a href="#">24075430</a>
KingFisher™ Deep-Well 96 Plate, V-bottom	<a href="#">95040450</a>
KingFisher™ 96 tip comb for DW magnets	<a href="#">97002534</a>
<b>MagMAX™ Express-96 instrument and accessories</b>	
MagMAX™ Express-96 Deep Well Magnetic Particle Processor	Contact your local sales representative.
MagMAX™ Express-96 Deep Well Plates	
MagMAX™ Express-96 Deep Well Tip Combs	

## Materials for PCR detection and analysis

**Table 3** TaqMan™ 2019-nCoV Assay Kit v1 (Cat. No. A47532)

Contents	Dye	Amount <sup>[1]</sup>	Concentration	Storage
2019-nCoV (ORF1ab) (Tube 1)	FAM™ dye	75 µL	20X	-30°C to -10°C
2019-nCoV (S Protein) (Tube 2)	FAM™ dye	75 µL	20X	
2019-nCoV (N Protein) (Tube 3)	FAM™ dye	75 µL	20X	
RNase P Assay (Tube 4)	VIC™ dye	250 µL	20X	

<sup>[1]</sup> Sufficient for 50 × 25-µL reactions.

**Table 4** TaqMan™ 2019-nCoV Control Kit v1 (Cat. No. A47533)

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

<sup>[1]</sup> Sufficient for 50 positive-control reactions.

**Table 5** Other required materials

Item	Source
<b>Master Mix</b>	
Thermo Scientific™ RNA UltraSense™ One-Step qPCR Master Mix: <ul style="list-style-type: none"> <li>• 250 µL RNA UltraSense™ Enzyme Mix</li> <li>• 1 mL RNA UltraSense™ 5X Reaction Mix</li> <li>• 300 µL 20X Bovine Serum Albumin (BSA)</li> <li>• 1 mL 50-mM Magnesium Sulfate (MgSO<sub>4</sub>)</li> <li>• 100 µL ROX™ Reference Dye</li> </ul>	11732927
Total RNA Control (Human)	4307281, or equivalent
<b>Real-Time PCR Instrument (one of the following)</b>	
Applied Biosystems™ QuantStudio™ 5 Food Safety Real-Time PCR System	A36328 or A36320
Applied Biosystems™ 7500 Fast Food Safety Real-Time PCR System	A30299 or A30304

Table 5 Other required materials (continued)

Item	Source
<b>Real-Time PCR Instrument Software (one of the following)<sup>[1]</sup></b>	
Applied Biosystems™ QuantStudio™ Design and Analysis Software v1.5.1 or later	Included with QuantStudio™ 5 Real-Time PCR System
Applied Biosystems™ 7500 Software SDS v.1.4.2 or later	Included with 7500 Real-Time PCR Instrument
<b>Equipment</b>	
Laboratory freezers -30°C to -10°C and ≤-70°C	MLS
Centrifuge, with a rotor that accommodates deepwell microplates	
Microcentrifuge	
Laboratory mixer, vortex or equivalent	
Single and multichannel adjustable pipettors (1 µL to 1,000 µL)	
Cold block (96-well or 384-well) or ice	
<b>Tubes, plates, and other consumables</b>	
MicroAmp™ Fast Optical 96-Well Reaction Plate, 0.1 mL	<a href="#">4346907</a>
MicroAmp™ Optical Adhesive Film Kit	<a href="#">4313663</a>
MicroAmp™ Optical 8-Cap Strips	<a href="#">4323032</a>
MicroAmp™ Fast 8-Tube Strip, 0.1 mL	<a href="#">4358293</a>
Disposable gloves	MLS
Micropipette tips, aerosol resistant	
Swabs with synthetic tip and aluminium or plastic shaft, preferably large macrofoam tipped applicator	
<b>Reagents</b>	
Thermo Scientific™ Oxoid™ Phosphate Buffered Saline	<a href="#">BR0014G</a>
Ethanol, 95%	MLS
Isopropanol, 100%	
Nuclease-free water	

<sup>[1]</sup> Real-Time PCR Instrument Software is used to control the instrument and to collect instrument run data.

# Procedural guidelines

## Guidelines for sample collection and storage

### Sample collection:

- Pre-moisten swab (preferably a large macrofoam tipped applicator with plastic shaft) with Phosphate Buffered Saline (PBS).
- When sampling, apply pressure with the wet swab onto the surface, move in at least two different directions while rotating the swab stick. Avoid letting the swab dry completely.
- Place the swab back into the transport tube containing 2 mL of PBS.

---

**Note:** Viral Transport Media (VTM) can be used instead of PBS according to World Health Organization protocol on *Surface sampling of coronavirus disease (COVID-19): A practical "how to" protocol for health care and public health professionals*, Version: 1.1, February 2020 ([www.who.int](http://www.who.int)).

---

### Storage:

- Swabs should be placed at refrigeration temperature (2-8°C) no more than 15 minutes post-sampling.

---

**Note:** Analyze the swabs in the coming 24 hours as much as possible. If the swabs are not likely to be analyzed within 48 hours maximum, store the swabs preferably between -70°C and -80°C, and shipped on dry ice.

---

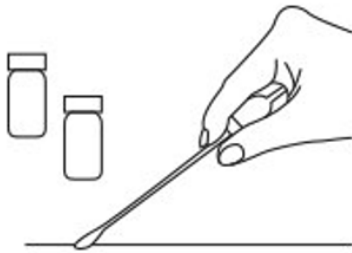
## Guidelines for the 2019-nCoV assay

- Use purified, non-degraded total nucleic acid that is free of RNase activity and RT-PCR inhibitors.
- Protect the assays and master mix from light.
- For each research sample, include the primers and probes for all three 2019-nCoV targets (FAM™ assay) in multiplex with the RNase P target (VIC™ assay).
- Before you begin, determine the number of required reactions. In addition to the nucleic acid research samples, include the following reactions for each 2019-nCoV assay:
  - One negative extraction control per plate
  - One 2019-nCoV Control v1 reaction per plate
  - One no-template control (NTC) per plate



## Workflow overview

### SARS-CoV-2 Detection on Food Packaging and Environmental Surfaces



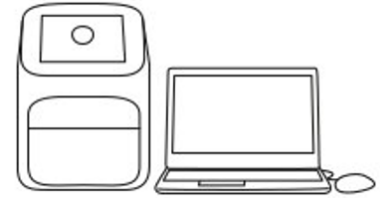
#### Sampling and Transport

Simple sample collection with Nylon swabs and Viral Transport Medium or Phosphate Buffered Saline



#### Sample Preparation

Options for automated medium to high-throughput and manual low-throughput



#### PCR and Data Analysis

Master mix and Real-Time PCR instrumentation with software designed to run environmental PCR range of products

## Prepare the samples and extract RNA using PrepSEQ™ Nucleic Acid Extraction Kit for Food and Environmental Testing

The KingFisher™ Flex-96 Deep-Well Magnetic Particle Processor or the MagMAX™ Express-96 Deep Well Magnetic Particle Processor can be used to extract RNA with the PrepSEQ™ Nucleic Acid Extraction Kit for Food and Environmental Testing.

- Ensure that the KingFisher™ Flex-96 Deep-Well Magnetic Particle Processor is set up with the KingFisher™ Flex -96 Deep-Well Heating Block.
- Ensure that the **PSNA\_Flex\_300ul** has been downloaded from the product page and loaded onto the instrument.

---

**Note:** Use **PSNA\_MagMAX\_300ul** if you are using MagMAX™ Express-96 Deep Well Magnetic Particle Processor.

---

- Add 74 mL 95% ethanol the Wash Buffer Concentrate bottle.
- Add 35 mL 100% isopropanol to an empty Binding Solution bottle.

1. Prepare and label the side of the processing plates according to the following table.

Plate ID	Plate type	Reagent	Volume per well
Wash Plate 1	KingFisher™ Deep-Well 96 Plates	Wash Buffer	300 µL
Wash Plate 2		Wash Buffer	300 µL
Elution Plate		Elution Buffer	100 µL
Tip Comb Plate	Place a KingFisher™ 96 tip comb for Deep Well magnets in a KingFisher™ 96 microplate		

2. Cover the plates with an adhesive film, then store at room temperature while setting up the sample plate (up to 1 hour).
3. Vortex the Magnetic Beads to ensure the bead mixture is homogenous.  
If a white precipitate has formed, first incubate the tube of Magnetic Particles at 37±1°C for approximately 10 minutes.

4. Prepare the Binding Bead Mix according to the following table.

Component	Volume per well <sup>[1]</sup>
Proteinase K Buffer	340 µL
Binding Solution	325 µL
Total Nucleic Acid Magnetic Beads	25 µL
Proteinase K	10 µL
<b>Total Binding Bead Mix</b>	<b>700 µL</b>

<sup>[1]</sup> Include 10% overage.

Prepare only the amount of Binding Bead Mix required for one day of use.

5. Invert the Binding Bead Mix five times gently to mix, then add 700 µL to each sample well and the Negative Extraction Control well in the Sample Plate.

---

**Note:** Remix the Binding Bead Mix by inversion frequently during pipetting to ensure even distribution of beads at all times. The Binding Bead Mix is viscous, so pipet slowly to ensure that the correct amount is added. DO NOT reuse pipette tips to add Binding Bead Mix to the samples, as the high viscosity will cause variations in the volumes added.

---

6. Add 1 µL of Total Human RNA Control to each sample well and to the Negative Extraction Control well.

7. Vigorously mix each sample tube containing the swab by vortexing for 30 seconds.

8. Add 300 µL of sample to each sample well and 300 µL of Nuclease-free Water (not DEPC-treated) to the Negative Extraction Control well in the Sample Plate.

9. Select the **PSNA\_Flex\_300ul** on the KingFisher™ Flex-96 Deep-Well Magnetic Particle Processor.

---

**Note:** Use **PSNA\_MagMAX\_300ul** if you are using MagMAX™ Express-96 Deep Well Magnetic Particle Processor.

---

10. Start the run, then load the prepared plates into position when prompted by the instrument.

11. After the run is complete (~40 minutes after start), immediately remove the Elution Plate from the instrument, then cover the plate with an adhesive film.

12. Place the Elution Plate on ice for immediate use in real-time RT-PCR.

---

**Note:** RNA can be stored at -70°C for long term storage (up to 1 year).

---

## Perform 1-Step RT-PCR

1. Set up the real-time PCR instrument as indicated in the following table

Stage	Step	Temperature	Time
Hold	Reverse transcription	50°C	15 minutes
Hold	Activation	95°C	2 minutes
Cycling (45 cycles)	Denaturation	95°C	3 seconds
	Anneal/Extension	60°C	30 seconds
Hold	(Optional) Melt curve	Refer to Real-Time instrument documentation.	

2. On ice, prepare a Reaction Mix for the number of reactions required plus 10% overage.

Component	Volume per reaction
RNA UltraSense™ Enzyme Mix	1 µL
RNA UltraSense™ 5X Reaction Mix	4 µL
2019-nCoV ORF1ab Assay (20X)	1 µL
2019-nCoV N Protein Assay (20X)	1 µL
2019-nCoV S Protein Assay (20X)	1 µL
RNase P Assay (20X)	1 µL
MgSO <sub>4</sub> (50 mM)	0.25 µL
<b>Total Reaction Mix Volume</b>	<b>9.25 µL</b>

3. For each reaction, combine the following components.

Component	Volume per reaction
Reaction Mix	9.25 µL
Template OR NEC OR 1 µL 2019-nCoV Control v1 + 9.75 µL Nuclease-free Water (not DEPC-Treated) OR NTC	10.75 µL
<b>Total Reaction Volume</b>	<b>20 µL</b>

4. Cap or seal the reaction vessels, and gently mix to make sure that all components are at the bottom of the amplification tube. Centrifuge briefly if needed.

---

**Note:** It is essential that no bubbles remain in the PCR reaction tubes. If bubbles remain after centrifugation, gently tap the reaction plate/tubes on the bench top and then briefly spin again at a higher speed until all bubbles are removed.

---

5. Place reactions in a preheated thermal cycler programmed as described above. Collect data and analyze results.

The following software should be used with each instrument.

Instrument	Software
7500 Fast Food Safety Real-Time PCR System	7500 Software SDS v.1.4.2 or later
QuantStudio™ 5 Food Safety Real-Time PCR System	QuantStudio™ Design and Analysis Software v1.5.1 or later

6. After cycling, hold the reaction at 4°C until further analysis.

## Analyze data

For more information about using a software, see the software user guide or help.

1. Open the data file (EDS) in the same software used to run the PCR.
2. Perform analysis using the following analysis settings.

Target	Baseline	Threshold
2019-nCoV assay (FAM™ dye)	Auto	Auto
RNaseP assay (IPC) (VIC™ dye)		

---

**IMPORTANT!** Analyze the run with no reference dye.

---

- For each plate, confirm that the control reactions for each 2019-nCoV assay perform as expected.

**Table 6** Expected results of the control reactions for each 2019-nCoV assay

Reaction	Expected $C_t$ value		About unexpected results
	2019-nCoV assay (FAM™ dye)	RNaseP assay (IPC) (VIC™ dye)	
NTC	Undetermined		If the NTC has an amplification curve that crosses the threshold (false positive), sample contamination may have occurred. Repeat the test with new reagents, following good RT-PCR practices.
Negative extraction control	Undetermined	$C_t < 40$	If a $C_t$ value is detected in the 2019-nCoV assay, then contamination may have occurred during extraction. Repeat the test with new reagents.
2019-nCoV Control v1	$C_t \leq 35$	$C_t < 40$	If the $C_t$ value is above the expected result, repeat the test with new reagents.

- Review all of the results for the 2019-nCoV assay to ensure that all positive results are detected. Evaluate the overall shape of the amplification curves. A sigmoidal amplification curve indicates true amplification.
- Export the results.

## Review results

- Classify the results for each individual assay, according to the  $C_t$  values in the table:

2019-nCoV assay (FAM™ dye)	RNaseP assay (IPC) (VIC™ dye)	2019-nCoV assay result
$C_t < 40$	Any value <sup>[1]</sup>	Positive. Confirm the result as described in Step 2.
$C_t = \text{Undetermined}$	$C_t < 40$	Negative.
$C_t = \text{Undetermined}$	$C_t = \text{Undetermined or } C_t = 40$	Invalid. Re-purify the nucleic acid from the sample, then repeat the test.

<sup>[1]</sup> An RNase P positive result is expected in most reactions. However, when a strong positive signal is detected in the 2019-nCoV assay the RNase P assay can occasionally give a negative result. These samples should be treated as positive if the amplification curve for the 2019-nCoV assay appears normal.

- To confirm a positive result, rerun the PCR reaction using the extracted sample RNA with each individual 2019-nCoV assay.

Component	Volume per reaction
RNA UltraSense™ Enzyme Mix	1 $\mu\text{L}$
RNA UltraSense™ 5X Reaction Mix	4 $\mu\text{L}$

*(continued)*

Component	Volume per reaction
2019-nCoV Assay	1 µL
RNase P Assay (20X)	1 µL
MgSO <sub>4</sub> (50 mM)	0.25 µL
Sample (Elution for RNA Extraction)	12.75 µL
<b>Total Reaction Volume</b>	<b>20 µL</b>

**Note:** This will result in three duplex PCR reactions being set up per sample being confirmed. Each reaction contains the primers and probes for one 2019-nCoV target (either ORF1ab, N protein or S protein) and the primers and probes for the RNase P target.

- For each test sample, interpret the results using the table. It is recommended that each lab perform accuracy testing with appropriate samples to establish guidelines for interpreting results.

2019-nCoV assay results	Interpretation of results
Any two of the three assays are positive.	SARS-CoV-2 RNA is present.
Any one of the assays is positive in two different samples collected from the same site.	SARS-CoV-2 RNA is present.
All three of the assays are negative.	SARS-CoV-2 RNA is not present.

# Troubleshooting

Observation	Possible cause	Recommended action
Inhibition of downstream PCR, indicated by non-detection of IPC reaction	Magnetic Particles were in the Elution Plate.	Avoid disturbing the Magnetic Particles during transfer of eluted RNA to the lyophilized assay.  Avoid transfer of Magnetic Particles using one of the following methods (optional): <ul style="list-style-type: none"> <li>Place the Elution Plate on the 96-Well Magnetic Ring Stand during transfer of eluted DNA sample to the lyophilized assay.</li> <li>Spin the plate at maximum speed in a plate centrifuge for the equivalent of approximately 4,000 × g for approximately 30 seconds, to pellet the Magnetic Particles to the bottom of the plate.</li> </ul>
	Elution Plate contains other inhibitory substances.	Dilute the eluted RNA 1:5 with Nuclease-free Water to dilute PCR inhibitors, and repeat the assay. If PCR remains inhibited, repeat the sample preparation.
In positive control wells, no target-specific signal is detected.	Positive control was omitted (pipetting error).	Repeat the assay. Make sure to pipet the positive control into all positive control wells.
In positive control wells or unknown sample wells, no IPC signal is detected, but target-specific signal is detected.	A high copy number of target RNA exists in samples, resulting in preferential amplification of the target-specific RNA.	No action is required. The result is considered positive if the target-specific amplification curve appears normal.
In negative extraction control wells, a target-specific signal is detected.	Carryover contamination occurred.	<ol style="list-style-type: none"> <li>Repeat the assay using fresh aliquots of all reagents and clean pipetting equipment.</li> <li>If the negative extraction control continues to show contamination, repeat the assay using a new kit.</li> <li>If the negative extraction control continues to show contamination, contact Technical Support.</li> </ol>
Amplification curve appears abnormal.	Auto baseline settings cannot account for large deviations in C <sub>T</sub> values between samples.	Manually adjust the baseline settings for the sample under investigation.
	Inhibitory substances carried over from RNA extraction.	Dilute the eluted RNA 1:5 with Nuclease-free Water to dilute PCR inhibitors, and repeat the assay. If PCR remains inhibited, repeat the sample preparation.



Observation	Possible cause	Recommended action
Amplification curve appears abnormal. <i>(continued)</i>	Primers and probes have degraded.	Repeat the PCR assay using fresh reagents.



# Supplemental information








## Good laboratory practices for PCR

To avoid amplicon contamination of samples, follow these guidelines when preparing or handling samples for PCR amplification:

- Wear clean gloves and a clean lab coat (not previously worn while handling amplified products or used during sample preparation).
- Change gloves whenever you suspect that they are contaminated.
- Maintain separate areas and dedicated equipment and supplies for:
  - Sample preparation and reaction setup.
  - Amplification and analysis of products.
- Do not bring amplified products into the reaction setup area.
- Open and close all sample tubes carefully. Avoid splashing or spraying samples.
- Keep reactions and components capped as much as possible.
- Use a positive-displacement pipettor or aerosol-resistant barrier pipette tips.
- Do not open reaction tubes after PCR.
- Do not autoclave reaction tubes after PCR.
- Clean lab benches and equipment periodically with 10% bleach solution or DNAZap™ Solutions (Cat. No. AM9890).

For additional information, refer to EN ISO 22174:2005 or [www.thermofisher.com/us/en/home/life-science/pcr/real-time-learning-center/real-time-pcr-basics.html](http://www.thermofisher.com/us/en/home/life-science/pcr/real-time-learning-center/real-time-pcr-basics.html).

## Symbol definitions

Symbol	Definition
	BATCH CODE
	CATALOG NUMBER
	CONTAINS SUFFICIENT FOR <n> TESTS
	CONSULT INSTRUCTIONS FOR USE
	MANUFACTURER
	UPPER AND LOWER TEMPERATURE LIMIT (storage temperature)
	USE BY



# Safety



**WARNING! GENERAL SAFETY.** Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
- Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, and so on). To obtain SDSs, see the “Documentation and Support” section in this document.

## Chemical safety



**WARNING! GENERAL CHEMICAL HANDLING.** To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below. Consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the "Documentation and Support" section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with sufficient ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if needed) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.



**WARNING! HAZARDOUS WASTE (from instruments).** Waste produced by the instrument is potentially hazardous. Follow the guidelines noted in the preceding General Chemical Handling warning.



**WARNING! 4L Reagent and Waste Bottle Safety.** Four-liter reagent and waste bottles can crack and leak. Each 4-liter bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position.

## Biological hazard safety



**WARNING! Potential Biohazard.** Depending on the samples used on this instrument, the surface may be considered a biohazard. Use appropriate decontamination methods when working with biohazards.



**WARNING! BIOHAZARD.** Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Conduct all work in properly equipped facilities with the appropriate safety equipment (for example, physical containment devices). Safety equipment can also include items for personal protection, such as gloves, coats, gowns, shoe covers, boots, respirators, face shields, safety glasses, or goggles. Individuals should be trained according to applicable regulatory and company/ institution requirements before working with potentially biohazardous materials. Follow all applicable local, state/provincial, and/or national regulations. The following references provide general guidelines when handling biological samples in laboratory environment.

- U.S. Department of Health and Human Services, *Biosafety in Microbiological and Biomedical Laboratories (BMBL)*, 5th Edition, HHS Publication No. (CDC) 21-1112, Revised December 2009; found at:  
<https://www.cdc.gov/labs/pdf/CDC-BiosafetymicrobiologicalBiomedicalLaboratories-2009-P.pdf>
- World Health Organization, *Laboratory Biosafety Manual*, 3rd Edition, WHO/CDS/CSR/LYO/2004.11; found at:  
[www.who.int/csr/resources/publications/biosafety/Biosafety7.pdf](http://www.who.int/csr/resources/publications/biosafety/Biosafety7.pdf)



# Documentation and support

## Food Safety support

Website: [thermoscientific.com/foodmicro](https://thermoscientific.com/foodmicro) or [thermofisher.com/foodsafety](https://thermofisher.com/foodsafety)

Support email:

- Europe, Middle East, Africa: [microbiology.techsupport.uk@thermofisher.com](mailto:microbiology.techsupport.uk@thermofisher.com)
- North America: [microbiology@thermofisher.com](mailto:microbiology@thermofisher.com)

Phone: Visit [thermofisher.com/support](https://thermofisher.com/support), select the link for phone support, and select the appropriate country from the dropdown menu.

## Related documentation

Document	Publication Number
<i>TaqMan™ 2019-nCoV Control Kit v1 Product Information Sheet</i>	MAN0019097
<i>QuantStudio™ 3 and 5 Real-Time PCR Systems Installation, Use, and Maintenance Guide</i>	MAN0010407
<i>Applied Biosystems™ 7300/7500/7500 Fast Real-Time PCR System Installation and Maintenance Guide</i>	4378657
<i>Applied Biosystems™ 7500/7500 Fast Real-Time PCR System: Maintenance Guide</i>	4387777

## Reference

World Health Organization protocol on *Surface sampling of coronavirus disease (COVID-19): A practical "how to" protocol for health care and public health professionals*, Version: 1.1, February 2020 ([www.who.int](http://www.who.int))

