

# Workflow and performance of the TaqPath COVID-19 RNase P Combo Kit 2.0

Multiplex real-time RT-PCR test for the qualitative detection of nucleic acids from SARS-CoV-2 and RNase P internal control

## Introduction

The Applied Biosystems™ TaqPath™ COVID-19 RNase P Combo Kit 2.0 contains assays and controls for a real-time reverse transcription polymerase chain reaction (RT-PCR) test intended for the qualitative detection of nucleic acids from SARS-CoV-2 in nasopharyngeal swabs and anterior or mid-turbinate nasal swabs from individuals suspected of COVID-19 by their health care provider, as well as individuals without symptoms.

- Human RNase P gene serves as an endogenous control
- The multiplex RT-PCR testing solution can detect the RNA from the SARS-CoV-2 virus and RNase P in a single reaction well
- Storage temperature of  $-20^{\circ}\text{C}$  for all components



Figure 1. Components of the TaqPath COVID-19 RNase P Combo Kit 2.0 (Cat. No. A51333).

## Workflow and turnaround time

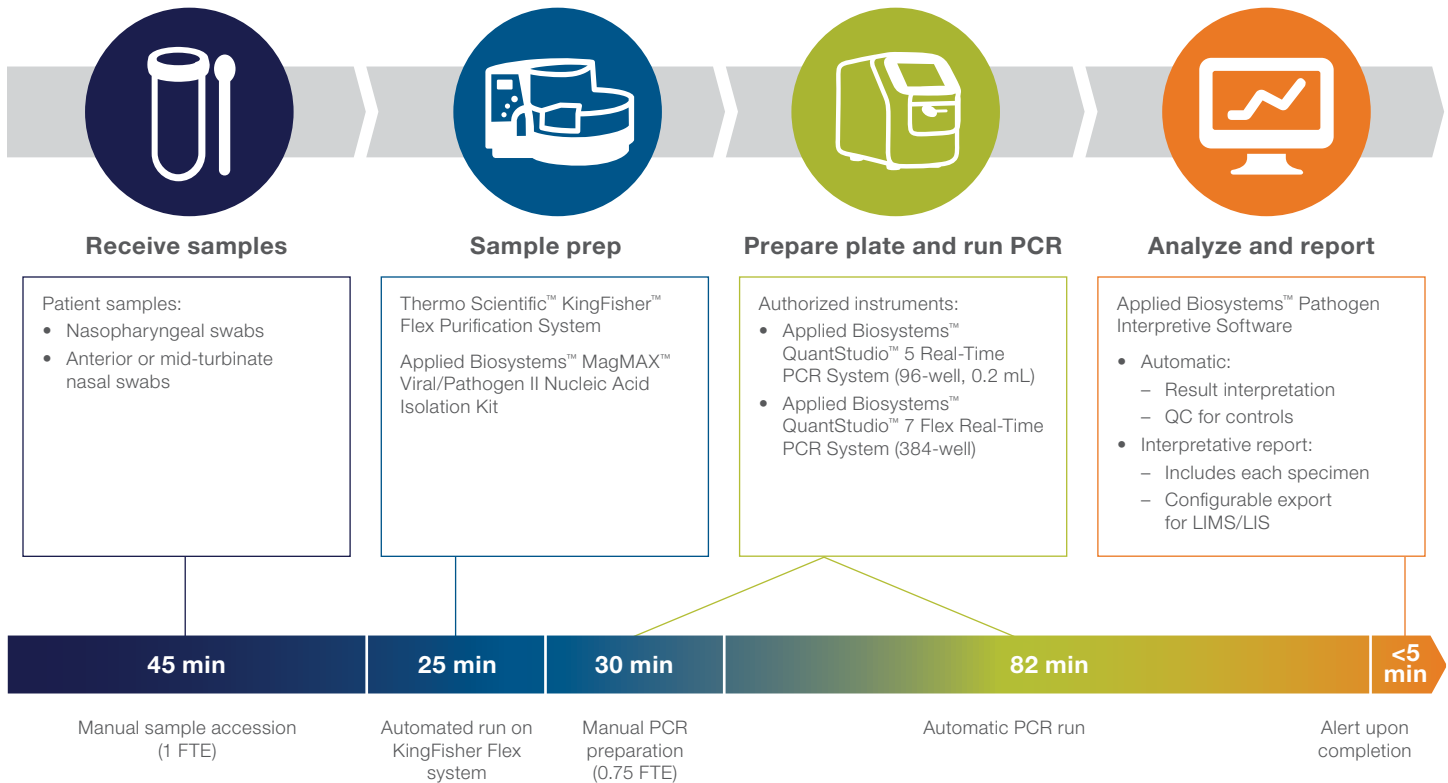


Figure 2. Workflow and turnaround time of the TaqPath COVID-19 RNase P Combo Kit.

The advanced assay design of the TaqPath COVID-19 RNase P Combo Kit compensates for current and future SARS-CoV-2 mutations:

- Unique fluorescence channel for each genomic region—*orf1a*, *orf1b*, and N genes (Figure 3)
- Redundancy with multiple targets (amplicons) per genomic region
- 8 targets spanning 3 genomic regions compensate for emerging mutations
- Excludes the S gene, which has a high risk of mutation
- Human RNase P gene serves as an endogenous specimen control

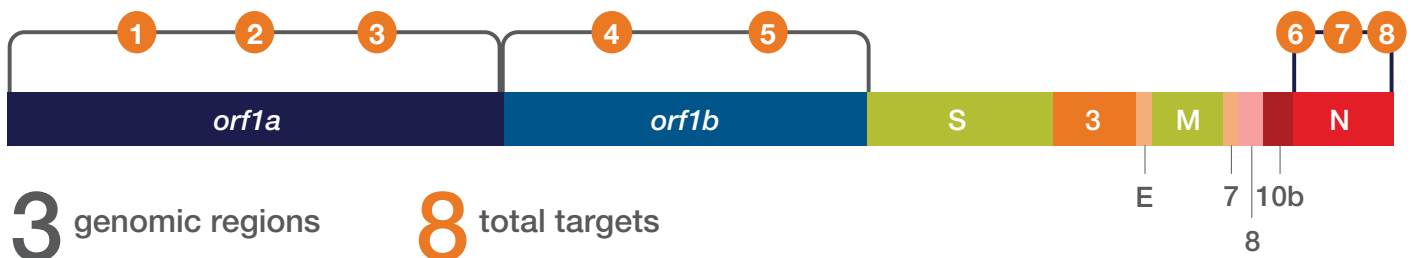


Figure 3. Schematic overview of the multi-target assay design of the TaqPath COVID-19 RNase P Combo Kit.

## Performance

### Limit of detection (LOD)

The LOD study established the lowest SARS-CoV-2 viral concentrations (genomic copy equivalents or GCE/mL) that can be detected at least 95% of the time using either anterior NS or NP swabs spiked with a gamma-irradiated SARS-CoV-2 isolate. The LODs for both specimen types were confirmed using 20 replicates each (Table 1).

### Reactivity (inclusivity)

*In silico* analysis was performed using 1,802,689 complete SARS-CoV-2 genomes in the GISAID and NCBI databases (June 9, 2021), including the variants of concern in Table 2.

All (100%) of the SARS-CoV-2 sequences of human origin exhibited 100% homology to one or more primer and probe sets, and are thus predicted to be detected.

### Interfering substances

The impact of potential interfering substances was tested by adding 10 potential interfering substances to pooled negative NP swab specimens spiked with gamma-irradiated SARS-CoV-2 at 3x the LOD (225 GCE/mL) and comparing them to a control with no interferents. Each interferent was tested in triplicate. No false-negative or false-positive results were observed for any interferents at the concentrations listed, on both the QuantStudio 5 and QuantStudio 7 Flex systems (Table 3).

**Table 1. LODs of Applied Biosystems™ QuantStudio™ PCR systems for both NS and NP swabs spiked with gamma-irradiated SARS-CoV-2.**

Real-time PCR instrument	Strain/isolate	LOD
QuantStudio 5 system (96-well, 0.2 mL)	USA-WA1/2020	75 GCE/mL NS
QuantStudio 7 Flex system (384-well)		75 GCE/mL NP swabs

**Table 2. *In silico* BLAST alignment to complete SARS-CoV-2 genomes.**

SARS-CoV-2 genomes	Number of genomes	Alignment
Human SARS-CoV-2	1,802,689	100%
Alpha variant	819,801	
Beta variant	20,057	
Gamma variant	25,068	
Delta variant	29,149	

**Table 3. Testing of interfering substances.**

Interfering substance	Concentration	Agreement
Mucin: bovine submaxillary gland, type I-S	0.1 mg/mL	100%
Blood (human)	1% v/v	100%
Afrin™ nasal spray	10% v/v	100%
Nasal corticosteroid—fluticasone propionate	5 µg/mL	100%
NeilMed™ Nasogel™ nasal gel	1% w/v	100%
bioAllers™ homeopathic allergy relief medicine	10% v/v	100%
Dorithricin™ throat lozenge, oral anesthetic and analgesic	1% w/v	100%
Oseltamivir phosphate	33 µg/mL	100%
Antibiotic, nasal ointment—pseudomonic acid	5 µg/mL	100%
Antibacterial, systemic—tobramycin	0.6 mg/mL	100%

## Cross-reactivity

### *In vitro*

Cross-reactivity testing was performed using the 31 organisms listed in Table 4. No false-positive SARS-CoV-2 calls were obtained with any organism tested.

### *In silico*

BLAST analysis was used to assess potential cross-reactivity between the TaqPath COVID-19 RNase P Combo Kit 2.0 primer/probe sequences and 55 organisms. With the exception of SARS-coronavirus, no bacterial, viral, or fungal microbe sequence aligned with  $\geq 80\%$  identity to more than one primer/probe set.

**Note:** The majority of the 285 SARS-coronavirus isolates shared  $\geq 80\%$  identity with more than one primer/probe set for the N gene and *orf1b*.

## Clinical evaluation

A clinical study was carried out to evaluate the performance of the TaqPath COVID-19 RNase P Combo Kit 2.0 using 120 archived nasopharyngeal swab and anterior nasal swab specimens.

The following specimens were tested:

- 60 samples that were positive for SARS-CoV-2 (54 nasopharyngeal swabs and 6 nasal swabs)
- 60 samples that were negative for SARS-CoV-2 (45 nasopharyngeal swabs and 15 nasal swabs)

The samples were tested using the TaqPath COVID-19 RNase P Combo Kit 2.0 and an FDA EUA-authorized comparator test. Positive percent agreement (PPA), negative percent agreement (NPA), and confidence intervals (CI) were calculated relative to the comparator test, summarized in Table 5.

**Table 4. Organisms used for *in vitro* cross-reactivity testing.**

Human coronavirus 229E	Influenza A	<i>Streptococcus pneumoniae</i>
Human coronavirus OC43	Influenza B	<i>Streptococcus pyogenes</i>
Human coronavirus HKU1	Enterovirus	<i>Bordetella pertussis</i>
Human coronavirus NL63	Respiratory syncytial virus	<i>Mycoplasma pneumoniae</i>
SARS-coronavirus*	Rhinovirus	<i>Pseudomonas aeruginosa</i>
MERS-coronavirus	Epstein-Barr virus	<i>Staphylococcus epidermidis</i>
Adenovirus	<i>Chlamydomphila pneumoniae</i>	<i>Streptococcus salivarius</i>
Human metapneumovirus	<i>Haemophilus influenzae</i>	<i>Candida albicans</i>
Parainfluenza 1	<i>Legionella pneumophila</i>	<i>Pneumocystis carinii</i>
Parainfluenza 2	<i>Mycobacterium tuberculosis</i>	Pooled human nasal wash
Parainfluenza 3		
Parainfluenza 4		

\* An inconclusive result with the SARS-coronavirus was obtained due to amplification of the N gene. Note: SARS-coronavirus is not a common respiratory pathogen and has not been in circulation since the 2003 outbreak.

**Table 5. Summary of clinical evaluation of the TaqPath COVID-19 RNase P Combo Kit 2.0 for nasal and nasopharyngeal sample types, compared to an FDA EUA-Authorized assay.**

Instrument	PPA	95% CI	NPA	95% CI
QuantStudio 5 system (96-well, 0.2 mL)	96.7%	88.5–99.6%	95.0%	86.1–99.0%
QuantStudio 7 Flex system (384-well)	95.0%	86.1–99.0%	96.7%	88.5–99.6%

## Conclusions

The TaqPath COVID-19 RNase P Combo Kit 2.0 is an ideal choice for COVID-19 testing using nasopharyngeal or anterior or mid-turbinate nasal swabs.

- Human sample confirmation
  - The RNase P gene serves as an endogenous control, helping to ensure sample integrity, quality, and extraction
- Innovative, multi-target primer/probe design
  - Eight targets across 3 regions (*orf1a*, *orf1b*, N gene) compensate for emerging SARS-CoV-2 mutations, providing confidence in results now and into the future
- Applied Biosystems Pathogen Interpretive Software
  - Helps decrease analysis and interpretation time, reducing the risk of user interpretation error
- Sensitive detection
  - Detects active infections in individuals suspected of having COVID-19 by their health care provider, as well as individuals without symptoms or other epidemiological reasons to suspect COVID-19
  - PCR-based test with excellent analytical sensitivity: 75 GCE/mL LOD
- Affordable and scalable
  - Increases testing throughput and lab efficiency
- Storage temperature of –20°C for all components

## Ordering information

Product	Components	Quantity	Cat. No.
TaqPath COVID-19 RNase P Combo Kit 2.0	TaqPath COVID-19 RNase P RT-PCR Kit 2.0 (1 tube of multiplex assay, 10 tubes of RNase P control)	1,000 reactions	A51333
	TaqPath COVID-19 Plus Control (10 tubes)		
	TaqPath COVID-19 Control Dilution Buffer (1 box with 10 tubes)		
TaqPath 1-Step Multiplex Master Mix (No ROX)	—	1 x 10 mL	A28523

Please refer to the Instructions for Use for more details and for the supported protocol. Deviations from the protocol are not permitted.

Find out more at [thermofisher.com/covid19evolved](https://thermofisher.com/covid19evolved)

**ThermoFisher**  
SCIENTIFIC