

PrimePCR SARS-CoV-2 Multiple and Single Mutation Assays



- SARS-CoV-2 surveillance by RT-qPCR — profile the mutation fingerprint of positive samples
- Rapid time to results — <60 minute protocol for use by researchers
- Flexible assay formats — multiple and single mutation detection
- High sensitivity — RT-qPCR assays can detect mutant alleles in samples at low RNA levels
- Expanding menu — assay options can be updated as new mutations and variants arise

Identify currently relevant mutations in severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) to rapidly screen samples for multiple variants of interest or concern, or survey the prevalence of a specific mutation using reverse transcription quantitative PCR (RT-qPCR) in a research use only (RUO) format.

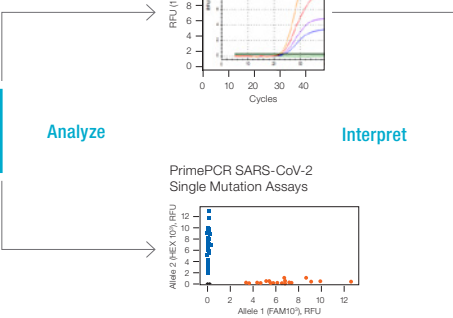
The PrimePCR SARS-CoV-2 Multiple Mutation Assays contain a primer set, up to six differentially labeled probes with homology to specific spike protein gene mutations, and an internal RNA integrity control. When used in combination with SARS-CoV-2 Standard and Variant Controls from Exact Diagnostics (a Bio-Rad Laboratories, Inc. company), the assays can be used by researchers to rapidly profile the molecular fingerprint of RNA from prior sample extractions. The SARS-CoV-2 Variant and SARS-CoV-2 Standard (wild-type) positive controls can be included to monitor reverse transcription, amplification, and detection steps. The high sensitivity of these assays can

be validated by researchers for use in assessing low-level samples with quantification cycle (Cq) values up to 40.

PrimePCR SARS-CoV-2 Single Mutation Assays target specific individual mutations in the SARS-CoV-2 genome that are believed to confer an evolutionary advantage to the virus. Assays contain a primer set plus two probes: one probe (labeled with HEX) matches the reference genome and the other probe (labeled with FAM) matches the mutation. The presence or absence of a specific mutation can be assessed in extracted SARS-CoV-2-positive samples using allelic discrimination. Each assay also contains a primer set and a probe as an internal control to assess the RNA integrity of the sample.

As part of an ongoing surveillance program, Bio-Rad can add new assays as novel mutations are acquired by the virus. Visit [bio-rad.com/SC2Variants](https://www.bio-rad.com/SC2Variants) to view the complete list of assays available.

Assay Workflows



Assay 1

Primer → ATTO 647 Cy5.5 HEX FAM Texas Red ← Primer

Amino acid change: K417N L452R MPC E484K/Q N501Y

SARS-CoV-2 Variant	K417N	L452R	MPC	E484K	E484Q	N501Y
Alpha (B.1.1.7)			*	*	*	*
Beta (B.1.351)	*	*	*	*	*	*
Delta (B.1.617.2)		*	*	*	*	*
Epsilon (B.1.427/B.1.429)		*	*	*	*	*
Gamma (P.1)		*	*	*	*	*
Kappa (B.1.617.1/ B.1.617.3)		*	*	*	*	*

Assay 2

Primer → ATTO 647 Cy5.5 HEX FAM Texas Red ← Primer

Amino acid change: K417T L452R MPC T478K N501Y

SARS-CoV-2 Variant	K417T	L452R	MPC	T478K	N501Y
Alpha (B.1.1.7)			*	*	*
Beta (B.1.351)	*	*	*	*	*
Delta (B.1.617.2)		*	*	*	*
Epsilon (B.1.427/B.1.429)		*	*	*	*
Gamma (P.1)	*	*	*	*	*
Kappa (B.1.617.1/ B.1.617.3)		*	*	*	*



Feature	PrimePCR SARS-CoV-2 Single Mutation Assay	PrimePCR SARS-CoV-2 Multiple Mutation Assay
Assay format (RUO)	Primer set with reference and mutant allele probes	Primer set with up to 6 probes
Mutations detected	1	Up to 5
Fluorophores	FAM HEX Texas Red	FAM ATTO 647 HEX Cy5.5 Texas Red
Internal RNA integrity control	Primer set and a probe to <i>RPP40</i>	Probe to a conserved region in spike gene between the primers (multiplexed control [MPC])
Protocol time	<60 minutes	
Product sizes available	200 reactions 1,000 reactions 2,500 reactions	

PrimePCR SARS-CoV-2 Multiple Mutation Assay Design

Mutation profiling can be used to assess the probability that a sample contains a SARS-CoV-2 variant of interest or concern. PrimePCR SARS-CoV-2 Multiple Mutation Assays include probes with homology to the sequences that encode the following key spike protein mutations:

Assay 1

SARS-CoV-2 Variant	K417N	L452R	E484K	E484Q	N501Y	MPC
Alpha (B.1.1.7)					•	•
Beta (B.1.351)	•		•		•	•
Delta (B.1.617.2)		•				•
Epsilon (B.1.427/B.1.429)		•				•
Gamma (P.1)			•		•	•
Kappa (B.1.617.1/ B.1.617.3)		•		•		•

MPC, multiplexed control.

Assay 2

SARS-CoV-2 Variant	K417T	L452R	T478K	N501Y	MPC
Alpha (B.1.1.7)				•	•
Beta (B.1.351)				•	•
Delta (B.1.617.2)		•	•		•
Epsilon (B.1.427/B.1.429)		•			•
Gamma (P.1)	•			•	•
Kappa (B.1.617.1/ B.1.617.3)		•			•

MPC, multiplexed control.

Related Products

Catalog # Description

Single Mutation Assays

12016270 PrimePCR SARS-CoV-2 Variant Assay, 200 reactions
 12016287 PrimePCR SARS-CoV-2 Variant Assay 1,000 reactions
 12016286 PrimePCR SARS-CoV-2 Variant Assay, 2,500 reactions

Multiple Mutation Assays

12016711 PrimePCR SARS-CoV-2 Variant Multi Kit, 200 reactions
 12016731 PrimePCR SARS-CoV-2 Variant Multi Kit, 1,000 reactions
 12016689 PrimePCR SARS-CoV-2 Variant Multi Kit, 2,500 reactions

Exact Diagnostics SARS-CoV-2 Reference Control (RUO)

COV019 SARS-CoV-2 Standard

Exact Diagnostics SARS-CoV-2 Variant Mutation Controls (RUO)

COVA SARS-CoV-2 S Gene Alpha Variant Control (B.1.1.7)
 COVB SARS-CoV-2 S Gene Beta Variant Control (B.1.351)
 COVE SARS-CoV-2 S Gene Epsilon Variant Control (B.1.427 and B.1.429)
 COVG SARS-CoV-2 S Gene Gamma Variant Control (P.1)

Exact Diagnostics SARS-CoV-2 Negative Control (RUO)

COV000 SARS-CoV-2 Negative

Instruments

12011319 CFX Opus 96 Real-Time PCR System
 12011452 CFX Opus 384 Real-Time PCR System
 1855195 CFX96 Touch Real-Time PCR System
 1855485 CFX384 Touch Real-Time PCR System
 1845097-IVD, CFX96 Dx Real-Time PCR System
 1841000-IVD

Reagents and PCR Plastics

12010176 Reliance One-Step Multiplex RT-qPCR Supermix, 1 ml, 200 x 20 µl reactions
 HSP9955 or equivalent* Hard-Shell 96-Well PCR Plates, low profile, thin wall, skirted, white shell/white well, barcoded, pkg of 50
 HSP3805 or equivalent* Hard-Shell 384-Well PCR Plates, low profile, thin wall, skirted, clear shell/white well, pkg of 50
 MSB1001 Microseal 'B' PCR Plate Sealing Film, adhesive, optical, pkg of 100

* Refer to the Hard-Shell PCR Plates Brochure (bulletin 5496) for other 96- and 384-well colored shell/white well PCR plates.



Visit bio-rad.com/SC2Variants for more information.



Bio-Rad PCR reagents and analytical instruments are manufactured under an ISO 13485:2016 certified Quality Management System and are quality control tested to ensure consistent product performance you can trust.

BIO-RAD, HARD-SHELL, and MICROSEAL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

The PrimePCR SARS-CoV-2 Multiple and Single Mutation Assays are for research use only. Not for use in diagnostic procedures.



Bio-Rad Laboratories, Inc.

Life Science Group

Website bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399
 Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23
 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050
 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670
 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23
 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23
 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

