



# Reliance SARS-CoV-2 Assays

## One-Step Multiplex RT-PCR

### Sensitive

- Established, gold-standard 1-step RT-PCR workflow for reliable detection of viral RNA

### Multiplexed

- Proven performance for the detection of multiple viral targets in a single reaction

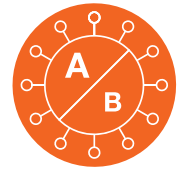
### Compatible

- Validated with most common RT-PCR systems

### Convenient

- Includes molecular controls

Introducing two molecular in vitro diagnostic tests from Bio-Rad for the detection of SARS-CoV-2: the Reliance SARS CoV-2 RT-PCR Assay Kit, for the qualitative detection of nucleic acid from SARS-CoV-2, and the Reliance SARS-CoV-2/FluA/FluB RT-PCR Assay Kit, for the detection and differentiation of SARS-CoV-2, influenza A, and influenza B. These kits are authorized for use with clinical samples under an Emergency Use Authorization (EUA) granted by the U.S. Food and Drug Administration (FDA).



### These kits include all reagents required to perform an RT-PCR test:

- Independently developed positive and negative molecular controls from Exact Diagnostics
- Oligonucleotide primers and probes, using sequences designed by the U.S. Centers for Disease Control and Prevention (CDC)
- Reliance One-Step Multiplex RT-qPCR Supermix, optimized for sensitive amplification of multiple targets in a single reaction

Both assays target the SARS-CoV-2 nucleocapsid (N) gene, which has higher viral specificity and may have lower mutation risk than the spike protein gene. Variant analysis confirms that these SARS-CoV-2 assays are not affected by current circulating escape variants (UK, South Africa, Brazil, and California variants), as of February 2021.

### The streamlined qPCR workflow can be finished in less than two hours\*



\* Excludes sample extraction time.

**Features of the Reliance SARS-CoV-2 RT-PCR Assay Kits**

Features	Reliance SARS-CoV-2 RT-PCR Assay Kit	Reliance SARS-CoV-2/FluA/FluB RT-PCR Assay Kit
Intended use*	Qualitative detection of nucleic acid from SARS-CoV-2 in specimens	Simultaneous, qualitative detection and differentiation of SARS-CoV-2, influenza A, and influenza B viral nucleic acid in specimens
Sensitivity (limit of detection)	125–500 copies/ml	SARS-CoV-2: 125–250 copies/ml Influenza A: 501–1,002 copies/ml Influenza B: 862 copies/ml
Targets detected	SARS-CoV-2 nucleocapsid (N) gene (N1 and N2 regions) Human RNaseP (internal control)	SARS-CoV-2: nucleocapsid (N) gene Influenza A: matrix (M1) gene Influenza B: nonstructural 2 (NS2) gene Human RNaseP (internal control)
Instrument compatibility	Bio-Rad CFX Opus 96 Real-Time PCR System Bio-Rad CFX Opus 384 Real-Time PCR System Bio-Rad CFX96 Touch Real-Time PCR System Bio-Rad CFX384 Touch Real-Time PCR System Bio-Rad CFX96 Dx System Applied Biosystems 7500 Fast Real-Time PCR System	Bio-Rad CFX Opus 96 Real-Time PCR System Bio-Rad CFX96 Touch Real-Time PCR System Bio-Rad CFX96 Dx System Applied Biosystems 7500 Fast Dx Real-Time PCR Instrument
Validated sample extraction methods	Thermo Fisher Scientific MagMAX Viral/Pathogen Nucleic Acid Isolation Kit QIAGEN QIAamp Viral RNA Mini Kit	QIAGEN QIAamp Viral RNA Mini Kit
Throughput	Up to 1,524 samples per day on a 384-well platform	Up to 372 samples per day on a 96-well platform

\* Refer to product instructions for use (IFU) for additional information.

**Ordering Information**

Catalog #	Description
12014115	Reliance SARS CoV-2 RT-PCR Assay Kit, 200 reactions
12015361	Reliance SARS CoV-2/FluA/FluB RT-PCR Assay Kit, 200 reactions

**Related Products**

**Reagents**

012010176	Reliance One-Step Multiplex RT-qPCR Supermix, 200 x 20 µl reactions, 1 ml
COV000	Exact Diagnostics SARS-CoV-2 Negative, 5 x 0.3 ml
COV019	Exact Diagnostics SARS-CoV-2 Standard, 5 x 0.3 ml
COVFLU	Exact Diagnostics SARS-CoV-2, Flu, RSV Run Control Positive, 5 x 1 ml
COVFLUNEG	Exact Diagnostics SARS-CoV-2, Flu, RSV Run Control Negative, 5 x 1 ml

**Instruments**

12011319	CFX Opus 96 Real-Time PCR System, 96-well, 5-color plus FRET, network-connected real-time PCR detection system, includes CFX Opus 96 base unit, cables
12011452	CFX Opus 384 Real-Time PCR System, 384-well, 5-color plus FRET, network-connected real-time PCR detection system, includes CFX Opus 384 base unit, cables
1855195	CFX96 Touch Real-Time PCR System, modular thermal cycler platform, includes C1000 Touch Thermal Cycler Chassis, CFX96 Optical Reaction Module, cables
1855485	CFX384 Touch Real-Time PCR System, modular thermal cycler platform, includes C1000 Touch Thermal Cycler Chassis, CFX384 Optical Reaction Module, cables
1845097-IVD, 1841000-IVD	CFX96 Dx Real-Time PCR System, CFX96 Dx optical reaction module and C1000 Dx Thermal Cycler

**PCR Plastics**

HSP9955*	Hard-Shell 96-Well PCR Plates, low profile, thin wall, skirted, white/white, pkg of 50
HSP3805*	Hard-Shell 384-Well PCR Plates, low profile, thin wall, skirted, clear/white, pkg of 50
MSB1001	Microseal 'B' PCR Plate Sealing Film, adhesive, optical, pkg of 100

\* Or equivalent. Refer to the Hard-Shell PCR Plate Brochure (Bio-Rad bulletin 5496) for other 96- and 384-well colored shell/white well PCR plates.

Visit [bio-rad.com/RelianceCOVIDKits](http://bio-rad.com/RelianceCOVIDKits) 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.

- These products have not been FDA cleared or approved but have been authorized for emergency use by the FDA under an EUA
- Reliance SARS CoV-2 RT-PCR Assay has been authorized only for the detection of nucleic acid from SARS-CoV-2, and not for any other viruses or pathogens
- Reliance SARS CoV-2/FluA/FluB RT-PCR Assay has been authorized only for the simultaneous qualitative detection and differentiation of nucleic acids from SARS-CoV-2, influenza A virus, influenza B virus and not for any other viruses or pathogens
- The emergency use of these products is only authorized for the duration of the declaration that circumstances justifying the authorization of emergency use of in vitro diagnostics for detection and/or diagnosis of COVID-19 under Section 564(b)(1) of the Act, 21 U.S.C. § 360bbb-3 (b)(1), unless the declaration is terminated, or authorization is revoked sooner

BIO-RAD, HARD-SHELL, and MICROSEAL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

Hard-Shell Plates are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 7,347,977; 6,340,589; and 6,528,302.

All trademarks used herein are the copyright of their respective owner.



**Bio-Rad  
Laboratories, Inc.**

---

Life Science  
Group

**Website** [bio-rad.com](http://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

