



Reliance SARS-CoV-2 Assay

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 the Reliance SARS-CoV-2 RT-PCR Assay Kit, a molecular in vitro diagnostic test for the qualitative detection of nucleic acid from SARS-CoV-2. The kit is authorized for use with clinical samples under an Emergency Use Authorization (EUA) granted by the U.S. Food and Drug Administration (FDA).



This kit includes 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

The assay targets the SARS-CoV-2 nucleocapsid (*N*) gene, which has higher viral specificity and may have lower mutation risk than the spike protein gene.

Coming soon: 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 (currently under FDA review for EUA).

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 Kit

Features	Reliance SARS-CoV-2 RT-PCR Assay Kit
Intended use*	Qualitative detection of nucleic acid from SARS-CoV-2 in upper respiratory tract specimens
Sensitivity (limit of detection)	125–500 copies/ml
Targets detected	SARS-CoV-2 nucleocapsid (N) gene (two targets) Human RNaseP (RP; 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
Validated sample extraction methods	Thermo Fisher Scientific MagMAX Viral/Pathogen Nucleic Acid Isolation Kit QIAGEN QIAamp Viral RNA Mini Kit
Throughput	Up to 1,524 samples per day on a 384-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

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

Instruments

12011319	CFX Opus 96 Real-Time PCR System , 96-well, 5-color plus fluorescence resonance energy transfer (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 shell/white well, pkg of 50
HSP3805*	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

* 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 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.

- This product has not been FDA cleared or approved but has been authorized for emergency use by FDA under an EUA for use by authorized laboratories
- The product has been authorized only for the detection of nucleic acid from SARS-CoV-2, not for any other viruses or pathogens
- The emergency use of this product is only authorized for the duration of the declaration that circumstances exist 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 Federal Food, Drug, and Cosmetic 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 **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

