

FACT SHEET FOR HEALTHCARE PROVIDERS

Bio-Rad Reliance SARS-CoV-2 RT-PCR Assay Kit – Bio-Rad Laboratories, Inc.

August 18, 2020

**Coronavirus
Disease 2019
(COVID-19)**

This Fact Sheet informs you of the significant known and potential risks and benefits of the emergency use of the Bio-Rad Reliance SARS-CoV-2 RT-PCR Assay Kit.

The Bio-Rad Reliance SARS-CoV-2 RT-PCR Assay Kit is authorized for use on upper respiratory specimens collected from individuals suspected of COVID-19 by their healthcare provider.

All patients whose specimens are tested with this assay will receive the Fact Sheet for Patients: Bio-Rad Reliance SARS-CoV-2 RT-PCR Assay Kit.

What are the symptoms of COVID-19?

Many patients with confirmed COVID-19 have developed fever and/or symptoms of acute respiratory illness (for example, cough, difficulty breathing). However, limited information is currently available to characterize the full spectrum of clinical illness associated with COVID-19. Based on what is known about the virus that causes COVID-19, signs and symptoms may appear any time from 2 to 14 days after exposure to the virus. Based on preliminary data, the median incubation period is approximately 5 days, but may range from 2 to 14 days.

Public health officials have identified cases of COVID-19 infection throughout the world, including the United States, which may pose risks for public health. Please check the Centers for Disease Control and Prevention (CDC) **Coronavirus (COVID-19) web page** for the most up-to-date information.

What do I need to know about COVID-19 testing?

Current information about COVID-19 for healthcare providers is available at CDC's web page. *Information for Healthcare Professionals* (see the links provided in "Where can I go for updates and more information?" section).

This test is to be performed using only respiratory specimens collected from individuals suspected of COVID-19 by their healthcare provider.

- The Bio-Rad Reliance SARS-CoV-2 RT-PCR Assay Kit can be used to test nasopharyngeal swabs, oropharyngeal swabs, nasal swabs, anterior nasal swabs, mid-turbinate nasal swabs, nasal aspirates, and nasal washes.
- The Bio-Rad Reliance SARS-CoV-2 RT-PCR Assay Kit should be ordered for the detection of COVID-19 in individuals suspected of having COVID-19 by their healthcare provider.
- The Bio-Rad Reliance SARS-CoV-2 RT-PCR Assay Kit is only authorized for use in laboratories certified under the Clinical Laboratory Improvement Amendments of 1988 (CLIA), 42 U.S.C. §263a, to perform high complexity tests.

Specimens should be collected following appropriate infection control precautions. Current guidance for COVID-19 infection control precautions are available at the CDC's website (see the links provided in "Where can I go for updates and more information?" section).

Use appropriate personal protective equipment when collecting and handling specimens from individuals suspected of having COVID-19 as outlined in the CDC *Interim Laboratory Biosafety Guidelines for Handling and Processing Specimens Associated with Coronavirus Disease 2019 (COVID-19)*. For additional information, refer to CDC *Interim Guidelines for Collecting, Handling, and Testing Clinical Specimens from Persons Under Investigation (PUIs) for Coronavirus Disease 2019 (COVID-19)* (see links provided in "Where can I go for updates and more information?" section).

What does it mean if the specimen tests positive for the virus that causes COVID-19?

Report adverse events, including problems with test performance or results, to MedWatch by submitting the online FDA Form 3500 (<https://www.accessdata.fda.gov/scripts/medwatch/index.cfm?action=reporting.home>) or by calling **1-800-FDA-1088**

FACT SHEET FOR HEALTHCARE PROVIDERS

Bio-Rad Reliance SARS-CoV-2 RT-PCR Assay Kit – Bio-Rad Laboratories, Inc.

August 18, 2020

Coronavirus
Disease 2019
(COVID-19)

A positive test result for COVID-19 indicates that RNA from SARS-CoV-2 was detected, and the patient is infected with the virus and presumed to be contagious. Laboratory test results should always be considered in the context of clinical observations and epidemiological data in making a final diagnosis and patient management decisions. Patient management should follow current CDC guidelines.

The Bio-Rad Reliance SARS-CoV-2 RT-PCR Assay Kit has been designed to minimize the likelihood of false positive test results. However, in the event of a false positive result, risks to patients could include the following: a recommendation for isolation of the patient, monitoring of household or other close contacts for symptoms, patient isolation that might limit contact with family or friends and may potentially increase contact with other patients with COVID-19, limits in the ability to work, the delayed diagnosis and treatment for the true infection causing the symptoms, unnecessary prescription of a treatment or therapy, or other unintended adverse effects.

All laboratories using this test must follow the standard testing and reporting guidelines according to their appropriate public health authorities.

What does it mean if the specimen tests negative for the virus that causes COVID-19?

A negative test result means that SARS-CoV-2 RNA was not present in the specimen above the limit of detection. However, a negative result does not rule out COVID-19 and should not be used as the sole basis for treatment or patient management decisions. A negative result does not exclude the possibility of COVID-19.

When diagnostic testing is negative, the possibility of a false negative result should be considered in the context of a patient's recent exposures and the presence of clinical signs and symptoms consistent with COVID-19. The possibility of a false negative result should especially be considered if the patient's recent exposures or clinical presentation indicate that COVID-19 is likely, and diagnostic tests for other causes of

illness (for example, other respiratory illness) are negative. If COVID-19 is still suspected based on exposure history together with other clinical findings, re-testing should be considered by healthcare providers in consultation with public health authorities.

Risks to a patient of a false-negative result include: delayed or lack of supportive treatment, lack of monitoring of infected individuals and their household or other close contacts for symptoms resulting in increased risk of spread of COVID-19 within the community, or other unintended adverse events.

The performance of this test was established based on the evaluation of a limited number of clinical specimens. The clinical performance has not been established in all circulating variants but is anticipated to be reflective of the prevalent variants in circulation at the time and location of the clinical evaluation. Performance at the time of testing may vary depending on the variants circulating including newly emerging strains of SARS-CoV-2 and their prevalence, which change over time.

What is an EUA?

The United States (U.S.) FDA has made this test available under an emergency access mechanism called an Emergency Use Authorization (EUA). The EUA for this test is supported by the Secretary of the Department of Health and Human Service's (HHS's) declaration that circumstances exist to justify the emergency use of in vitro diagnostics (IVD) for the detection and/or diagnosis of the virus that causes COVID-19.

An IVD made available under an EUA has not undergone the same type of review as an FDA-approved or cleared IVD. The FDA may issue an EUA when certain criteria are met, which includes that there are no adequate, approved, available alternatives, and based on the totality of scientific evidence available, it is reasonable to believe that this IVD may be effective in the detection of the virus that causes COVID-19.

The EUA for this test is in effect for the duration of the COVID-19 declaration justifying the emergency use of

Report adverse events, including problems with test performance or results, to MedWatch by submitting the online FDA Form 3500 (<https://www.accessdata.fda.gov/scripts/medwatch/index.cfm?action=reporting.home>) or by calling **1-800-FDA-1088**

FACT SHEET FOR HEALTHCARE PROVIDERS

Bio-Rad Reliance SARS-CoV-2 RT-PCR Assay Kit – Bio-Rad Laboratories, Inc.
August 18, 2020

Coronavirus
Disease 2019
(COVID-19)

IVD, unless it is terminated or revoked by the FDA (after which the test may no longer be used).

Where can I go for updates and more information?

CDC webpages:

General: [cdc.gov/COVID19](https://www.cdc.gov/COVID19)

Healthcare Professionals:

[cdc.gov/coronavirus/2019-nCoV/guidance-hcp.html](https://www.cdc.gov/coronavirus/2019-nCoV/guidance-hcp.html)

Information for Laboratories: [cdc.gov/coronavirus/2019-nCoV/guidance-laboratories.html](https://www.cdc.gov/coronavirus/2019-nCoV/guidance-laboratories.html)

Laboratory Biosafety: [cdc.gov/coronavirus/2019-nCoV/lab-biosafety-guidelines.html](https://www.cdc.gov/coronavirus/2019-nCoV/lab-biosafety-guidelines.html)

Isolation Precautions in Healthcare Settings:

[cdc.gov/coronavirus/2019-ncov/infection-control/control-recommendations.html](https://www.cdc.gov/coronavirus/2019-ncov/infection-control/control-recommendations.html)

Specimen Collection: [cdc.gov/coronavirus/2019-nCoV/guidelines-clinical-specimens.html](https://www.cdc.gov/coronavirus/2019-nCoV/guidelines-clinical-specimens.html)

Infection Control: [cdc.gov/coronavirus/2019-ncov/infection-control/index.html](https://www.cdc.gov/coronavirus/2019-ncov/infection-control/index.html)

FDA webpages:

General: [fda.gov/novelcoronavirus](https://www.fda.gov/novelcoronavirus)

EUAs: (For links to patient fact sheet and manufacturer's instructions)

[fda.gov/medical-devices/emergency-situations-medical-devices/emergency-use-authorizations](https://www.fda.gov/medical-devices/emergency-situations-medical-devices/emergency-use-authorizations)

Bio-Rad, Laboratories, Inc.:

4000 Alfred Nobel Drive
Hercules, CA USA 94547

Phone: 510-741-1000

Fax: 510-741-5800

Website: [bio-rad.com](https://www.bio-rad.com)

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.

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

Report adverse events, including problems with test performance or results, to MedWatch by submitting the online FDA Form 3500 (<https://www.accessdata.fda.gov/scripts/medwatch/index.cfm?action=reporting.home>) or by calling **1-800-FDA-1088**