



QIAstat-Dx Respiratory SARS-CoV-2 Panel

For quick and accurate detection of respiratory infections



Influenza-like illnesses aren't just caused by influenza viruses. There are many possible bacterial and viral causes that result in similar, overlapping symptoms. As a result, diagnosing respiratory infections using traditional methods can be slow and challenging. Imagine getting 23 results in about an hour. Powered by our easy workflow and reliable multiplex PCR technology, the QIAstat-Dx Respiratory SARS-CoV-2 Panel makes this possible.

- **Get comprehensive results fast**

- Detect and differentiate 23 pathogen targets in about an hour
- Support timely clinical decision-making and rapid therapeutic intervention

- **Set up your tests quickly and easily**

- All-in-one assay cartridges require less than one minute hands-on time
- Dry swab or just 300 μ L liquid transport medium with no precision pipetting required

- **Receive supportive information**

- Easily access Ct values and amplification curves for all detected pathogens



Comprehensive pathogen coverage with QIAstat-Dx

Viral

- SARS-CoV-2
- Influenza A
- Influenza A subtype H1N1/2009
- Influenza A subtype H1
- Influenza A subtype H3
- Influenza B
- Coronavirus 229E
- Coronavirus HKU1
- Coronavirus NL63
- Coronavirus OC43
- Parainfluenza virus 1
- Parainfluenza virus 2
- Parainfluenza virus 3
- Parainfluenza virus 4
- Adenovirus
- Respiratory syncytial virus A/B
- Human metapneumovirus A/B
- Bocavirus
- Rhinovirus/Enterovirus*

Bacterial

- *Mycoplasma pneumoniae*
- *Legionella pneumophila*
- *Bordetella pertussis*
- *Chlamydomphila pneumoniae*

*Enterovirus and Rhinovirus are both detected, but not differentiated, with the QIAstat-Dx Respiratory SARS-CoV-2 Panel.



QIAstat-Dx Respiratory SARS-CoV-2 Panel runs on both our modular **QIAstat-Dx Analyzer** and higher-throughput **QIAstat-Dx Rise**.

Ordering Information - Assay

Product	Contents	Cat. no.
QIAstat-Dx Respiratory SARS-CoV-2 Panel	Six individually packaged cartridges containing all reagents needed for sample preparation and multiplex RT-real time PCR plus internal control, including six transfer pipettes	691214



Confidently detect respiratory infections with QIAstat-Dx – learn more at [QIAGEN.com/QIAstat-Dx](https://www.qiagen.com/QIAstat-Dx)

The QIAstat-Dx Analyzer, QIAstat-Dx Rise and QIAstat-Dx Respiratory SARS-CoV-2 Panel are intended for in vitro diagnostic use.

Trademarks: QIAGEN®, Sample to Insight®, QIAstat-Dx® (QIAGEN Group). PROM-15948-002 1127825 07/2022 © 2022 QIAGEN, all rights reserved.

Ordering
Technical Support
Website

www.qiagen.com/shop
www.support.qiagen.com
www.qiagen.com