

AMPIPROBE® SARS-CoV-2 Assay Kit (RUO)

Detection of SARS-CoV-2 RNA by RT-qPCR

Our AMPIPROBE® SARS-CoV-2 Kit (RUO) is a multiplexed assay designed for the detection of SARS-CoV-2 virus. It contains two primer/probe sets specific to different SARS-CoV-2 genomic regions and primers/probes for internal control, human RNase P, which is also present in the negative control. This kit is intended to be used in combination with general RNA extraction kits, including AMPIXTRACT™ SARS-CoV-2 Extraction Kit (RUO) and generic qPCR instruments. This assay can also be used in manual settings or with open automated platforms such as GENFLEX® platform (please contact us for more information) or any other open platforms available in the market for the liquid handling process of RNA extraction and PCR set up.

Ordering Information

[Order Online »](#)

ENZ-GEN230-0100	96Reactions
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

- **Reliable** – Accurate detection of SARS-CoV-2 RNA using our highly optimized AMPIPROBE® technology
- **Adaptable** – Adaptable to the high throughput GENFLEX™ platform (sample processing, amplification and detection) or any automated open platforms, and manual settings
- **Affordable** - Low-cost alternative to other methods of viral load detection
- **Validated:** Tested with several qPCR instruments such as QuantStudio™ 5 (Applied Biosystems), RotorGene (QIAGEN), q-Tower³ (Anlaytik Jena) and compatible with any qPCR instruments able to read 3 channels (Cy5, FAM, HEX)

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Coronavirus (COVID-19)
Application	qPCR
Application Notes	Detection of SARS-CoV-2 RNA by RT-qPCR
Contents	AMPIPROBE [®] RT-PCR Enzyme Mix AMPIPROBE [®] RT-PCR Buffer AMPIPROBE [®] SARS-CoV-2 Primer/Probe Mix Nuclease-free Water
Sensitivity	98.10%
Specificity	99.30%

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)