



Detection Kit for SARS-CoV-2

AMPIPROBE® SARS-CoV-2 Assay Kit

RUO

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 our 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.

GENFLEX® platform



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)

Affordable

Low-cost alternative to other detection methods of viral load detection

Validated

Tested with several qPCR instruments such as QuantStudio™ 5 (Applied Biosystems®), RotorGene® Q (QIAGEN®), q-Tower3 (Analytik Jena) and compatible with any qPCR instruments able to read 3 channels (Cy5, FAM, HEX)

AMPIPROBE® SARS-CoV-2 Assay Kit (RUO)

ENZ-GEN230-0100

COMPONENTS FOR 96 TESTS	QUANTITY	STORAGE
AMPIPROBE® RT-PCR Buffer	1.3 mL	-20°C
AMPIPROBE® SARS-CoV-2 Primer/Probe Mix	260 µL	-20°C
AMPIPROBE® RT-PCR Enzyme Mix	130 µL	-20°C
Nuclease-free Water	1 mL	-20°C

AMPIPROBE® SARS-CoV-2 Controls (RUO)

ENZ-GEN231-0004

The AMPIPROBE® SARS-CoV-2 Controls are designed to be used during the PCR process to ensure an accurate RNA detection and amplification.

COMPONENTS	QUANTITY	STORAGE
AMPIPROBE® SARS-CoV-2 Positive Control	100 µL	-80°C
AMPIPROBE® SARS-CoV-2 Negative Control	100 µL	-80°C
No Template Control (NTC)	100 µL	-80°C

RELATED PRODUCTS	
Product Name	Product Number
AMPIXTRACT® SARS-CoV-2 Extraction Kit (RUO)	ENZ-GEN225
SARS-CoV-2 IgG ELISA Kit (RUO)	ENZ-GEN190

Global Headquarters
ENZO LIFE SCIENCES, INC.
Ph: 800.942.0430
info-usa@enzolifesciences.com

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

**Belgium, The Netherlands
& Luxembourg**
Ph: +32 3 466 0420
info-be@enzolifesciences.com

France
Ph: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Ph: +49 7621 5500 526
info-de@enzolifesciences.com

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

For detailed product information, visit us online: www.enzolifesciences.com/coronavirus

For Research Use Only (RUO)

All product names, logos, and brands are the property of their respective owners.