

AMPIPROBE[®] SARS-CoV-2 Controls (RUO)

Controls for AMPIPROBE[®] SARS-CoV-2 Assay Kit

The AMPIPROBE[®] SARS-CoV-2 Controls (RUO) contains a positive RNA control that is specific to the SARS-CoV-2 genomic regions targeted by the assay as well as human RNase P, a negative RNA control that contains target to human RNase P, and a no-template control (NTC). This kit can be used in combination with AMPIPROBE[®] SARS-CoV-2 Assay Kit (RUO) and Controls (RUO).

Ordering Information

[Order Online »](#)

ENZ-GEN231-0004	1Kit
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- Includes positive, negative, and no template controls

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name Coronavirus (COVID-19)

Application qPCR

Application Notes SARS-CoV-2 RNA Controls for qPCR

Contents Positive control
Negative control
No template control

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)