

AMPIXTRACT™

SARS-CoV-2

Extraction Kit (RUO)

Extraction of SARS-CoV-2 RNA for qPCR

The AMPIXTRACT™ SARS-CoV-2 Extraction Kit (RUO) utilizes magnetic-particle technology for automated isolation and purification of nucleic acids from various biological specimens. This kit is intended to be used in combination AMPIPROBE® SARS-CoV-2 Assay kit (RUO) and Controls (RUO). The extraction can be performed manually or on Enzo's GENFLEX® platform (please contact us for more information) or on any extraction platform open to accept third party reagents.

Magnetic-particle technology enables purification of high-quality nucleic acids that are free of proteins, nucleases, and other impurities. The purified nucleic acids are ready for direct use in downstream applications, such as amplification and qualitative detection of microorganisms in biological specimens, enzymatic reactions or for storage for later use.

Please contact us if you are considering to use our extraction kit on alternative platforms.

Ordering Information

[Order Online »](#)

ENZ-GEN225-0100	96Reactions
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

- **Effective** - Isolate high-quality RNA using magnetic bead extraction and optimized extraction reagents.
- **Reliable** - Consistent RNA yield from a small amount of starting material
- **Validated** - Tested to work for downstream applications such as RT-qPCR

Handling & Storage

Use/Stability Upon receipt:(1) AMPIXTRACT™ Wash Reagent and AMPIXTRACT™ Lysis Reagent should be stored at room temperature.(2) AMPIXTRACT™ Elution Buffer, AMPIXTRACT™ Magnetic Beads, and AMPIXTRACT™ PK Dilution Buffer should be stored at 4°C.(3) AMPIXTRACT™ Stabilizer, AMPIXTRACT™ Carrier RNA, and AMPIXTRACT™ Proteinase should be -20°C.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Coronavirus (COVID-19)

Application qPCR

Application Notes Extraction of SARS-CoV-2 RNA for qPCR

Contents

- AMPIXTRACT™ Wash Reagent
- AMPIXTRACT™ Lysis Reagent
- AMPIXTRACT™ Elution Buffer
- AMPIXTRACT™ Magnetic Beads
- AMPIXTRACT™ PK Dilution Buffer
- AMPIXTRACT™ Stabilizer
- AMPIXTRACT™ Proteinase
- AMPIXTRACT™ Carrier RNA

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-
fr@enzolifesciences.com](mailto:info-fr@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)