

AMPIXTRACT® SARS-CoV-2 Extraction Kit

RUO

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 with AMPIPROBE® SARS-CoV-2 Assay and Controls kits (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 bead 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.

GENFLEX® platform



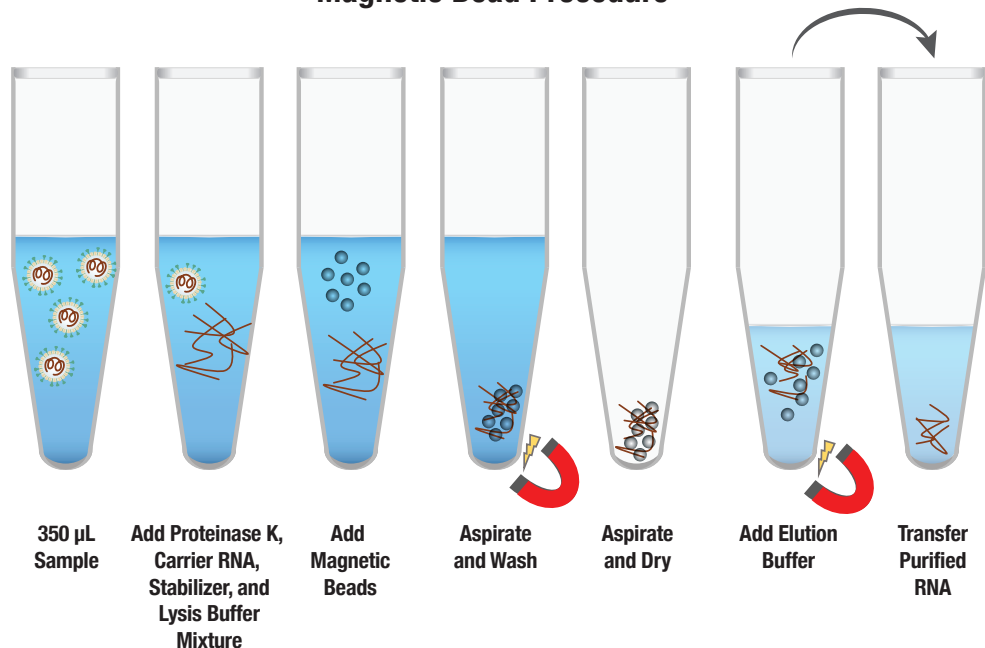
High Performance

- Magnetic bead separation technology with proprietary buffer system
- Isolation of high-quality RNA in < 1 hour
- > 99% recovery of extracted RNA from 350 µL sample

Easily Adaptable

- Manual or automated processing
- Compatible with common automated open platforms and RT-PCR kits
- Flexible for 1.5 mL tubes or 96-well plates

Magnetic Bead Procedure



AMPIXTRACT® SARS-CoV-2 Extraction Kit (RUO)

ENZ-GEN225-0100

COMPONENTS FOR 96 TESTS	VOLUME	STORAGE
Wash Buffer	1 x 50 mL	RT (20-25°C)
Lysis Buffer	1 x 80 mL	RT (20-25°C)
Elution Buffer	1 x 12 mL	2-8°C
Magnetic Beads	1 x 5 mL	2-8°C
PK Dilution Buffer	1 x 9 mL	2-8°C
Stabilizer	2 x 900 µL	-20°C
Proteinase	1 x 350 µL	-20°C
Carrier RNA	1 x 1.2 mL	-20°C

RELATED PRODUCTS	
Product Name	Product Number
AMPIPROBE® SARS-CoV-2 Controls (RUO)	ENZ-GEN231
AMPIPROBE® SARS-CoV-2 Assay Kit (RUO)	ENZ-GEN230

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.