



Extraction Kit for Viral & Pathogen DNA and RNA

AMPIXTRACT® Virus & Pathogen Extraction Kit

Over 99% recovery of nucleic acids in less than one hour.

The AMPIXTRACT® Virus & Pathogen Extraction Kit utilizes magnetic-particle technology for optimized isolation and purification of nucleic acids from various biological specimens. This kit is intended to be used for downstream nucleic acid detection applications such as RT-qPCR.

The extraction can be performed either manually or on any automated extraction platform open to accept third-party reagents.

Magnetic bead technology enables the 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 quantitative detection of microorganisms in biological specimens, enzymatic reactions, or for storage for later use.

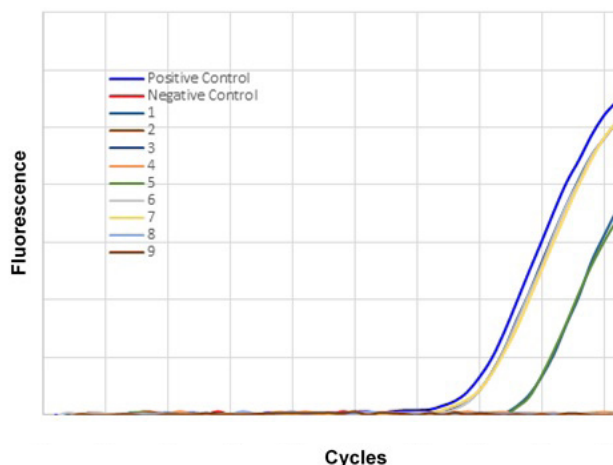
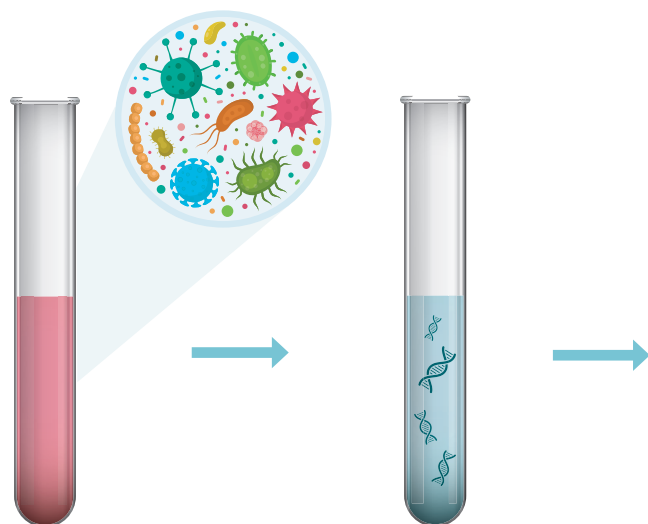
High Performance

- >99% recovery of extracted RNA in 1 hour
- Predictable and reproducible yields
- Magnetic bead separation technology with a proprietary buffer system

Easily Adaptable

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

Optimized virus and pathogen detection by qPCR



Extraction

AMPIXTRACT® Virus &
Pathogen Extraction Kit

Amplification

AMPIGENE® DNA polymerases

Detection

Dye & Probe based mixes



AMPIGENE® ladders

AMPIXTRACT® Virus & Pathogen Extraction Kit ENZ-GEN249

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	PRODUCT #	PRODUCT	PRODUCT #
AMPIGENE® qPCR 1-Step Green Kit Lo-ROX	ENZ-NUC108	CYGREEN® Nucleic Acid Dye	ENZ-GEN105
AMPIGENE® qPCR 1-Step Green Kit Hi-ROX	ENZ-NUC109	AMPIGENE® Taq Polymerase	ENZ-PRT100
AMPIGENE® qPCR 1-Step Probe Kit Lo-ROX	ENZ-NUC110	AMPIGENE® HS Taq DNA Polymerase	ENZ-PRT-101
AMPIGENE® qPCR 1-Step Probe Kit Hi-ROX	ENZ-NUC111	AMPIGENE® Taq Mix	ENZ-NUC100
AMPIGENE® qPCR 1-Step Probe No-ROX	ENZ-NUC112	AMPIGENE® HS Taq Mix	ENZ-NUC101
AMPIGENE® qPCR Green Mix Lo-ROX	ENZ-NUC103	AMPIGENE® dNTP Mix	ENZ-NUC102
AMPIGENE® qPCR Green Mix Hi-ROX	ENZ-NUC104	AMPIGENE® cDNA Synthesis Kit	ENZ-KIT106
AMPIGENE® qPCR Probe Mix Lo-ROX	ENZ-NUC105	AMPIGENE® 1-Step PT-PCR Kit	ENZ-KIT105
AMPIGENE® qPCR Probe Mix Hi-ROX	ENZ-NUC106		
AMPIGENE® qPCR Probe Mix No-ROX	ENZ-NUC107		

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 instrument compatibility and detailed product information visit us online: <https://www.enzolifesciences.com/pcr-qpcr>

Research Use Only. Not for Diagnostic Procedure.

FL012-Rev02282022