

Viral RNA+DNA Preparation Kit, L pack

Spin column based RNA+DNA purification from blood, serum, tissue, cell culture or swabs

Viral RNA+DNA Preparation Kit is designed for rapid and effective isolation of RNA and DNA from a variety of pathogen organisms such as virus or bacteria. Samples can be fresh or frozen plasma/blood (treated with anticoagulants excepted heparin), serum, other cell-free body fluids or pathogen-infected tissue. The kit allows high yield isolation of viral RNA/DNA from nasal or throat swabs. The Kit is specifically designed to isolate high-quality nucleic acids using low elution volumes and allowing sensitive downstream analysis including quantitative PCR and RT-PCR. The purified RNA/DNA is free of proteins and nucleases. Viral RNA+DNA Preparation Kit uses lysis buffer including chaotropic salts to inactivate RNases/DNases and advanced silica-gel membrane technology for fast purification of intact RNA/DNA. The preparation procedure is optimized to achieve reliable results within 30 min.

Ordering Information

[Order Online »](#)

JBS-PP-235L	250preparations
-------------	-----------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

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

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

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