

Viral RNA Extraction Kits

Isolate RNA from saliva or nasal/nasopharyngeal swabs

Isolate Viral RNA In Minutes

EpiGentek's **EpiQuik™ Viral RNA Isolation Kits** allow researchers to quickly prepare viral RNA directly from cell-free liquid specimens, specifically from saliva and nasal or nasopharyngeal swabs. The RNA purified with these kits can be used for a variety of routine applications.

EpiQuik™ Viral RNA Isolation Fast Kit

Cat #P-9107-050 (50 samples)

Rapid extraction of total viral RNA from saliva and nasal/nasopharyngeal swabs via spin columns

- Delivers high-quality RNA in 10 minutes
- Recover quantities of up to 5 µg of purified RNA

EpiMag™ Viral RNA Isolation Kit

Cat #P-9108-050 (50 samples)

Magnetic bead-based extraction of total viral RNA from saliva and nasopharyngeal swabs

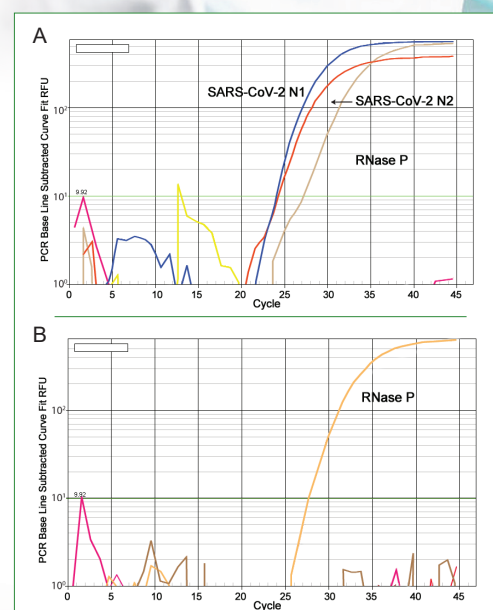
- Delivers high-quality RNA in 25 minutes
- Allows small size (>30 bases) and large size (<200 kb) to be isolated
- No need for centrifugation

EpiMag™ 96-Well Viral RNA Extraction Kit (High Throughput)

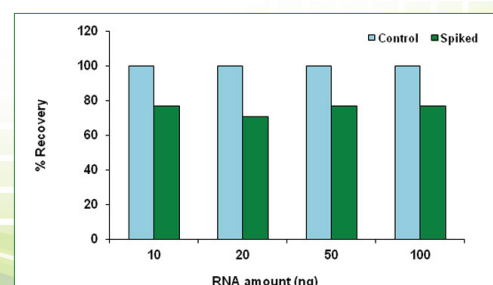
Cat #P-9109-96 (96 samples)

Magnetic bead-based high throughput extraction of total viral RNA from saliva and nasopharyngeal swabs

- Extract up to 96 samples simultaneously
- Suitable for robotic handling and automation setups
- Finish within 30 minutes
- No need for centrifugation



▲ Total RNA was isolated from nasal swab samples using the **EpiQuik Viral RNA Isolation Fast Kit** and used for RT-PCR analysis. A: Nasopharyngeal swab samples were spiked with SARS-CoV-2-N positive control; B: Nasopharyngeal swab samples were not spiked with SARS-CoV-2-N positive control. PCR was processed using primers and probes against SARS-CoV-2 N1/N2 and human RNase P gene (internal control).



▲ RNA at different concentrations was spiked into normal saline then isolated with the **EpiMag 96-Well Viral RNA Extraction Kit (High Throughput)** and fluorescently quantified (RNA/ssDNA).

Ordering Information

Product Name	Size	Cat. No.
EpiQuik Viral RNA Isolation Fast Kit	50 samples	P-9107-050
EpiMag Viral RNA Isolation Kit	50 samples	P-9108-050
EpiMag 96-Well Viral RNA Extraction Kit (High Throughput)	96 samples	P-9109-96



100% Satisfaction Guarantee

All EpiGentek products are covered by our 6-month money back guarantee

Learn more at:
www.epigentek.com/rna-isolation