

ViraBind™

Adenovirus Miniprep Kit

Purification of viruses via cesium chloride ultracentrifugation procedures can be tedious and time-consuming. ViraBind™ Adenovirus Purification Kits provide a much more efficient system for fast adenoviral purification with high yields. The ViraBind™ Adenovirus Miniprep Kit uses a special spin column to purify supernatant from a single T75 flask or 10cm dish per prep in about 30 minutes.

Ordering Information

[Order Online »](#)

| | |
|-------------|----------------|
| CBI-VPK-099 | 10preparations |
|-------------|----------------|

Manuals, SDS & CofA

[View Online »](#)

- Achieve the quality of CsCl procedures in <2 hours, without ultracentrifugation
- Recover >90% adenoviral yield

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