

QuickTiter™

Lentivirus Titer Kit (Lentivirus-Associated HIV p24), Trial Size

Measuring the HIV-1 p24 antigen is a long-established method for lentivirus quantitation. However, the traditional p24 ELISA detects both virus-associated p24 and free p24 generated by 293 cells during transient transfection. Free p24 can account for a substantial portion of total p24 in the supernatant. Therefore, the ELISA typically overestimates the quantity of lentivirus present. Our QuickTiter™ Lentivirus Titer Kit (Lentivirus-Associated HIV p24) substantially minimizes this problem. A proprietary technology separates the lentivirus from free p24 in solution prior to running the ELISA portion of the assay.

Ordering Information

[Order Online »](#)

CBI-VPK-107-T	32 assays
---------------	-----------

Manuals, SDS & CofA

[View Online »](#)

- Proprietary technology separates virus-associated p24 from free p24
- Minimizes overestimation of lentivirus titer common to traditional p24 ELISA kits
- Lentivirus quantitation on a standard microplate reader
- HIV-1 p24 Standard included

Handling & Storage

Use/Stability Upon receipt, aliquot and store the Recombinant HIV-1 p24 Standard at -20°C to avoid multiple freeze/thaw cycles. Store all other kit components at 4°C.

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

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](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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