

QuickTiter™ MuLV Core Antigen ELISA Kit

An anti-MuLV p30 monoclonal coating antibody is adsorbed onto a microtiter plate. MuLV core antigen (p30) present in the sample or standard binds to the antibodies adsorbed on the plate; an anti-MuLV p30 polyclonal antibody is added and binds to the antigen captured by the first antibody. Following incubation and wash steps, a HRP-conjugated secondary antibody is added and binds to the anti-MuLV p30 polyclonal. Unbound HRP-conjugated secondary antibody is removed during the wash steps, and substrate solution reactive with HRP is added to the wells. A colored product is formed in proportion to the amount of MuLV core antigen present in the sample. The reaction is terminated by addition of acid and absorbance is measured at 450 nm. A standard curve is prepared from recombinant MuLV p30 core antigen and sample concentration is then determined.

Ordering Information

[Order Online »](#)

CBI-VPK-156	96 assays
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

- Measures the MuLV core protein (p30) as low as 300 pg/mL
- Quantitation is performed on a standard microplate reader
- Recombinant MuLV p30 Standard included

Handling & Storage

Use/Stability Upon receipt, aliquot and store the Recombinant MuLV p30 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, ELISA

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)