

Ebola virus monoclonal antibody (FE25)

Ebola virus is an enveloped, negative-stranded RNA virus that causes disease in humans known as Ebola virus disease (EVD) or Ebola hemorrhagic fever (EHF). Symptoms typically begin with a fever, throat and muscle pains, and headaches approximately 2-21 days after contacting the virus. This is followed by nausea, vomiting and diarrhea, which is accompanied by decreased liver and kidney function, which can lead to bleeding. The virus is spread through contact with bodily fluids of an infected animal or person.

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

ENZ-ABS267-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Do not freeze.

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, WB

Application Notes Detects a band of ~30kDa by Western blot.

Clone FE25

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

Host Mouse

Immunogen Ebola virus (formalin inactivated, Zaire strain)

Isotype IgG2a

Purity Detail Protein G affinity purified.

Species Reactivity Ebola virus

Specificity Recognizes ebola virus in human samples.

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



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)