

# 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](#).

**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 <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



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)