

# Anthrax Lethal Factor monoclonal antibody (LF2)

Anthrax Lethal Factor (LF) is a zinc-dependent protease that cleaves mitogen-activated protein kinase kinase (MAPKK) and causes lysis of macrophages. LF, together with Protective Antigen, forms Lethal Toxin, which is the dominant virulence factor produced by *B. anthracis* and is the major cause of death in infected animals. Lethal toxin has been reported to trigger apoptosis in human melanoma cells.

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

## Ordering Information

[Order Online »](#)

ADI-905-621-100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** LF, ALF

**Application** ELISA, IP, WB

**Clone** LF2

**Formulation** Liquid. In PBS containing 0.01% sodium azide.  
Concentration 1mg/ml.

**Host** Mouse

**Immunogen** Anthrax lethal factor.

**Isotype** IgG1

**Purity Detail** Protein G affinity purified.

**Source** Purified from hybridoma tissue culture supernatant.

**Species Reactivity** Bacillus anthracis

**UniProt ID** P15917

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

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-be@enzolifesciences.com)

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