

Anthrax Protective Antigen monoclonal antibody (PA6)

The exotoxin produced by *B. anthracis* is composed of three proteins: a cell-receptor binding protein known as Protective Antigen (PA), and two enzymes, Edema Factor (EF) and Lethal Factor (LF). An 83 kDa form of PA binds to a cell surface receptor and, upon proteolytic activation to a 63 kDa fragment, heptamerizes to form a membrane channel that mediates the entry of three molecules of LF or EF into the cells. PA and LF form Lethal Toxin, which is the dominant virulence factor produced by *B. Anthracis* and is the major cause of death in infected animals.

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

Ordering Information

[Order Online »](#)

ADI-905-619-100	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

Application ELISA, IP, WB

Clone PA6

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

Host Mouse

Immunogen Anthrax protective antigen.

Isotype IgG1

Purity Detail Protein G affinity purified.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Bacillus anthracis

UniProt ID P13423

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-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@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)