

Aprotinin monoclonal antibody (21)

Aprotinin, also called bovine pancreatic trypsin inhibitor, is a single chain globular Kunitz-type serine proteinase inhibitor found in bovine pancreas, parotid gland and lung. It forms stable, enzymatically inactive complexes with trypsin, chymotrypsin, plasmin and kallikrein. Aprotinin has been used therapeutically as an antifibrinolytic agent under the name Trasylol (Bayer) to reduce blood loss during cardiopulmonary bypass surgery, but this use has been associated with an increased frequency of acute renal injury.

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

Ordering Information

[Order Online »](#)

BPD-ABS-037-21-1	1mg
BPD-ABS-037-21-04	400µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Bovine pancreatic trypsin inhibitor

Application ELISA, WB

Clone 21

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

GenBank ID 404172

Host Mouse

Immunogen Aprotinin from bovine lung.

Isotype IgG1κ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Bovine

Specificity Binds bovine aprotinin immobilized on a solid phase or free in solution.

UniProt ID P00974

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-
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)