

# Aprotinin monoclonal antibody (18)

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](#).

**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	18
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-  
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)

Last modified: May 29, 2024