

# Caspase-3 polyclonal antibody

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

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-210-806-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



**Figure:** Western blot analysis of human caspase-3 using PAb to caspase-3 (Prod. No. ALX-210-806) with (lane 1) and without (lane 2) blocking peptide.

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	WB
Application Notes	Detects bands of ~32kDa (procaspase-3) and ~20kDa (cleavage product) by Western blot.
Formulation	Liquid. In PBS, pH 7.2, containing 0.5% BSA, 5mM EDTA, 0.03% Proclin and 30% glycerol.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 163-175 of human caspase-3.
Purity Detail	Epitope-affinity purified.
Recommendation Dilutions/Conditions	Western Blot (0.5-4µg/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P42574
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)