

Caspase-1 polyclonal antibody

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

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-8040	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

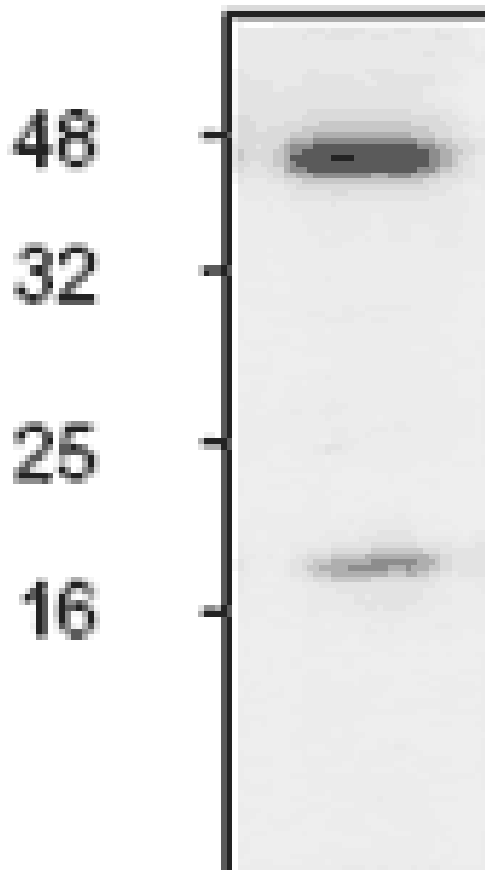
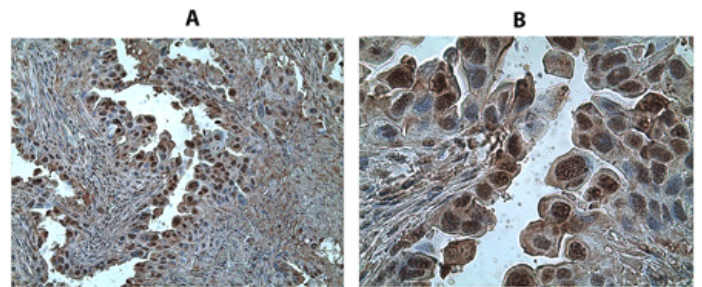


Figure: Western blot analysis of caspase-1 expression in rat liver tissue lysate using Caspase-1, pAb (Prod. No. [ALX-210-804](#)). Caspase-1 polyclonal antibody Western blot, Caspase-1 polyclonal antibody Western blot, yes

ALX-210-806, 210-806.gif, <http://static.enzolifesciences.com/fileadmin/files/image/210-806.gif>, **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.

ALX-210-806, 210-806.gif, **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.



Immunohistochemistry staining of human lung tumor tissue using Caspase-1, pAb (Prod. no. ALX-210-804) at a 1:50 dilution showed positive staining in both cytoplasm and nucleus of lung epithelial cells. Images were taken at magnification of 20 x (A) and 63 x (B).

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

Alternative Name	IL-1 β converting enzyme, Interleukin-1 β converting enzyme, ICE
Application	ICC, IHC, IP, WB
Application Notes	Detects bands of ~45kDa (procaspase-1) and ~20kDa (cleavage product) by Western blot.
Formulation	Liquid. In PBS, pH 7.2, containing 30% glycerol, 0.5% BSA, 0.5mM EDTA, and 0.03% Proclin.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 224-242 (CEPH224KTSDSTFLVFMSHGIREG242) of human caspase-1.
Purity Detail	Epitope-affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (10-20 μ g/ml)Immunohistochemistry (1:50)Immunoprecipitation (4-8 μ g/ml)Western Blot (0.5-4 μ g/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	P29466
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-
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)