

# Caspase-4 monoclonal antibody (4B9)

The caspases are a family of cysteine proteases that cleave after certain aspartate residues, and are primarily recognized as mediators of apoptosis. caspases are synthesized as inactive zymogens that can be cleaved to form active enzymes following the induction of apoptosis by stress or death receptors. Initiator caspases (e.g. caspase-8 and -10) are activated by dimerization of the zymogen on a dedicated adaptor protein. These activated initiator caspases in-turn cleave downstream effector or executioner caspases (e.g. caspase-3, -6, and -7) in a cascade-like manner, which cleave key cellular proteins that lead to the morphological changes associated with apoptotic cell death.

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

Citations: 9

[View Online »](#)

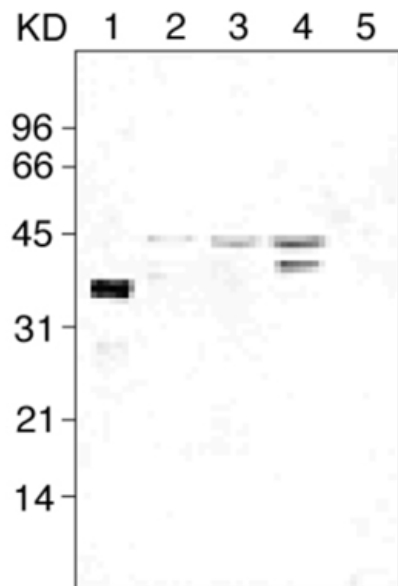
## Ordering Information

[Order Online »](#)

ADI-AAM-114-E	100µg
---------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of caspase-4 probed with Caspase-4, mAb (4B9). Lane 1: TXp20 (recombinant); Lane 2: U937 cell extract; Lane 3: HL60 cell extract; Lane 4: HUC-Fm cell extract; Lane 5: PC12 cell extract.

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Application      WB

Application Notes      Detects a band of ~45kDa by Western blot.

Clone      4B9

Formulation      Liquid. In PBS, pH 7.2, containing 50% glycerol.

GenBank ID      HSU28014

Host      Mouse

Immunogen      Recombinant human caspase-4 (N-terminal fragment).

Isotype      IgG1

Purity Detail      Protein A affinity purified.

**Recommendation**  
**Dilutions/Conditions**      Western Blot (1:1,000, ECL) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity      Human

UniProt ID      P49662

**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)