

# Caspase-9 monoclonal antibody (5B4)

Apoptosis is a major form of cell death characterized by several morphological features that include chromatin condensation and fragmentation, cell membrane blebbing, and formation of apoptotic bodies. These morphological changes occur via signalling pathways that lead to the recruitment and activation of caspases, a family of cysteine-containing, aspartate-specific proteases. Caspases exist as inactive proenzymes in cells, and are activated through their processing into two subunits in response to apoptotic stimulation. Activated caspases cleave a variety of important cellular proteins, other caspases, and Bcl-2 family members, leading to a commitment to cell death. Caspase-9, also known as ICE-LAP61, Mch-62 or Apaf-33 is a 45 kDa protein that is involved in one of the relatively well characterized caspase cascades. It is triggered by Cytochrome C release from the mitochondria, which promotes the activation of caspase-9 by forming a complex with Apaf-1 in the presence of dATP. Once activated, caspase-9 initiates a caspase cascade involving caspase-3, -6, and -7. The active form of caspase-9 is composed of a 32-35 kDa subunit and 10 kDa subunit. Caspase-9 is expressed in a variety of tissues.

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

Citations: 16

[View Online »](#)

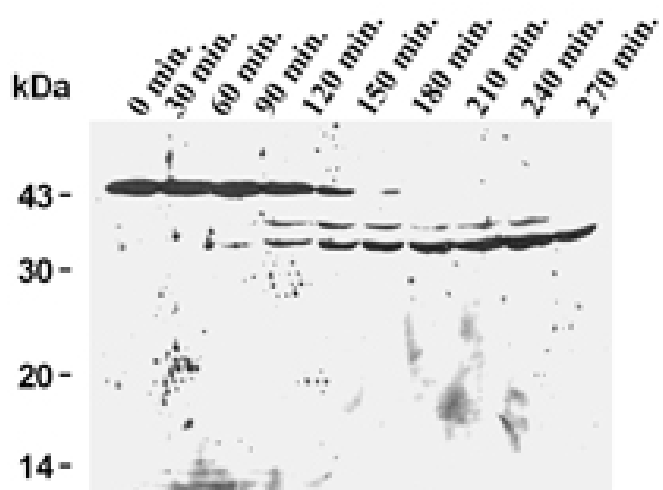
## Ordering Information

[Order Online »](#)

ADI-AAM-139-E	100µg
---------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Caspase-9 fragments expression in apoptosis induced Jurkat cells by Fas mAb, probed with Caspase-9 mAb (5B4).

## 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 ~46-50kDa by Western blot.
Clone	5B4
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol.
GenBank ID	AAC50640
Host	Mouse
Immunogen	Recombinant human caspase-9 (N-terminal fragment).
Isotype	IgG1
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Drosophila, Human, Mouse, Rat
UniProt ID	P55211
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



ENZO LIFE SCIENCES,  
Last modified **May 29, 2024**  
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-be@enzolifesciences.com](mailto:info-be@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)