

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: 17

[View Online »](#)

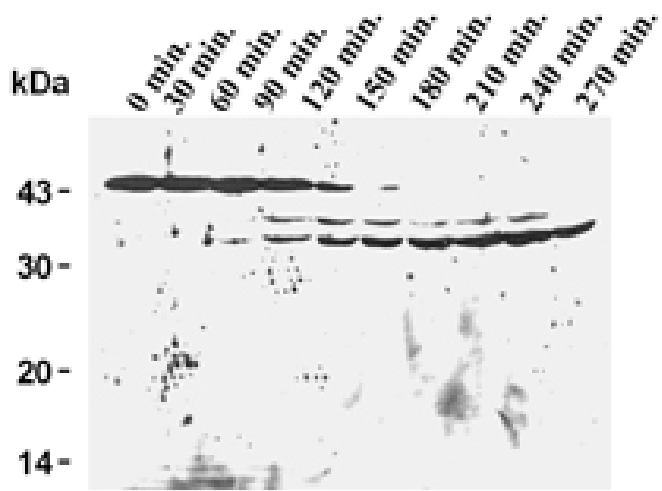
Ordering Information

[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 [Worry-Free Guarantee](#)



European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com