

Caspase-9 polyclonal antibody

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](#).

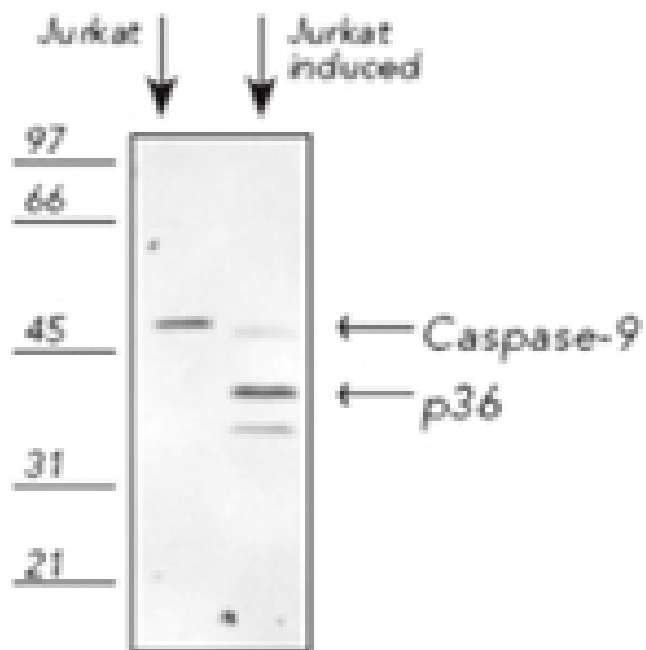
Ordering Information

[Order Online »](#)

ADI-AAP-149-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of untreated and apoptosisinduced Jurkat cell lysates, probed with Caspase-9 pAb.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application WB

Application Notes Detects bands of ~46-50kDa (proform) and ~35/37kDa, ~18kDa, ~10kDa (processed) by Western blot.

Formulation Liquid. In PBS containing 0.09% sodium azide.

GenBank ID AAC50640

Host Rabbit

Immunogen Recombinant human procaspase-9 (N-terminal fragment).

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:500, ECL) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Chicken, Dog, Guinea pig, Human, Monkey, Mouse, Rat, Sheep

UniProt ID P55211

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-
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)