

Caspase-3 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](#).

Citations: 6

[View Online »](#)

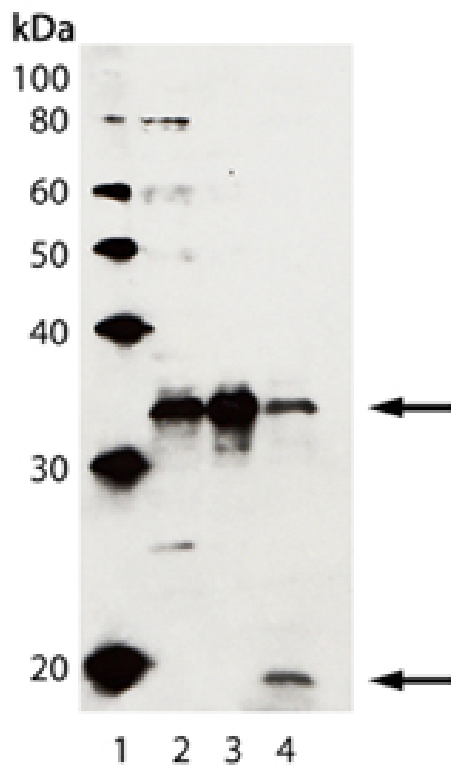
Ordering Information

[Order Online »](#)

ADI-AAS-103-D	50µl
ADI-AAS-103-F	200µl

Manuals, SDS & CofA

[View Online »](#)



Western Blot Analysis of Caspase-3, pAb (Prod. No. ADI-AAS-103): Lane 1:MW Marker; Lane 2: HeLa Cell Lysate (Prod. No. ADI-LYC-HL100); Lane 3: Jurkat Cell Lysate (Prod. No. ADI-LYC-JK100); Lane 4: Jurkat + Staurosporine

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 a band of ~36kDa by Western blot.

Formulation Liquid. Antiserum.

GenBank ID U26943

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of human caspase-3 . The sequence differs from rat caspase-3 by 1 aa and from mouse by 2 aa.

Recommendation Dilutions/Conditions Western Blot (ECL, 1:5,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Whole rabbit serum.

Species Reactivity Bovine, Chicken, Dog, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep, Xenopus

UniProt ID P42574

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



ENZO LIFE SCIENCES,
INC.
Phone: +1 940 948 9120
info@enzolifesciences.com
usa@enzolifesciences.com

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

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

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

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

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