

Caspase-10 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](#).

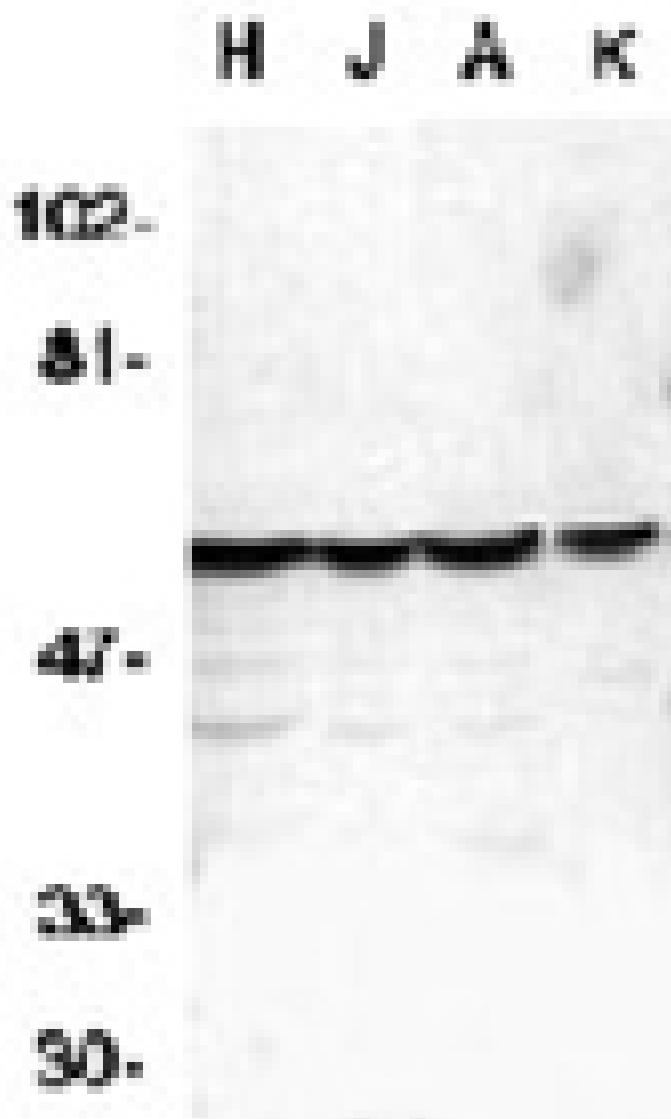
Ordering Information

[Order Online »](#)

ADI-AAP-110-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Caspase-10 in HeLa (H), Jurkat (J), A431 (A), K562 (K) whole cell lysates with Caspase-10 antibody at 1 μ g/ml.

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 ~59kDa by Western blot.

Formulation Liquid. In PBS containing 0.02% sodium azide.

GenBank ID U86214

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 505-521 of human Caspase-10.

Purity Detail DEAE purified.

Recommendation Dilutions/Conditions Western Blot (1.0µg/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human

UniProt ID Q92851

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

.

Last modified: May 29, 2024



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)