

Caspase-7 monoclonal antibody (10-1-62)

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.

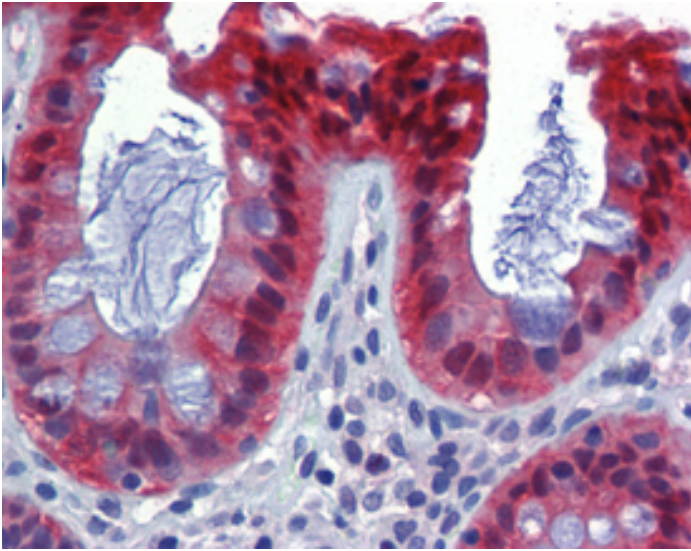
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

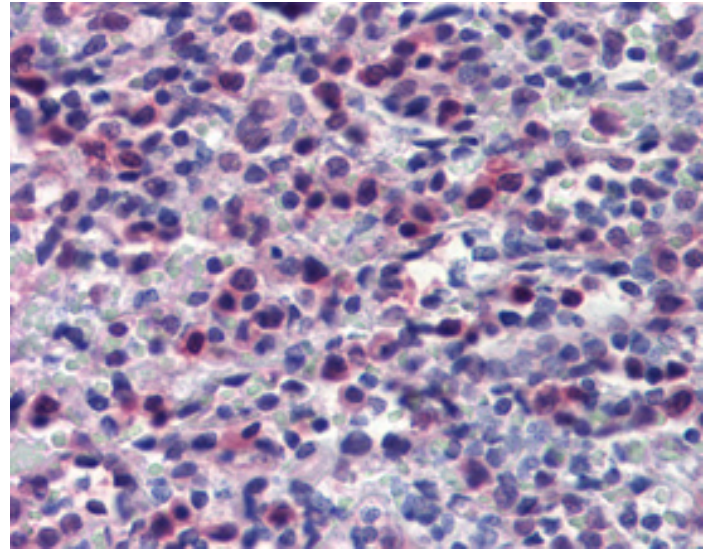
[View Online »](#)

Manuals, SDS & CofA

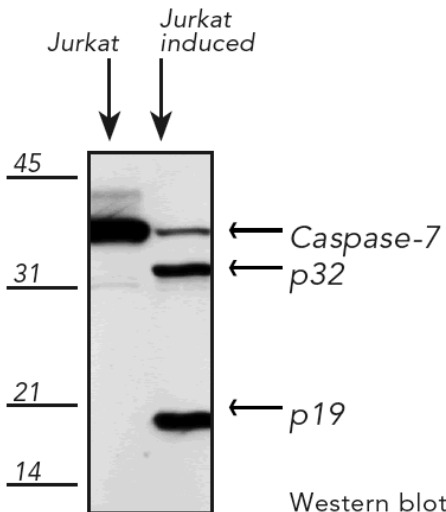
[View Online »](#)



IHC Analysis: Human colon tissue stained with Caspase-7, mAb (10-1-62) at 10µg/ml.



IHC Analysis: Human spleen tissue stained with Caspase-7, mAb (10-1-62) at 10µg/ml.



Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC (PS), WB

Application Notes Immunohistochemistry (paraffin sections)
Western Blot (1µg/ml, ECL)
Optimal conditions must be determined individually for each application.

Clone 10-1-62

Formulation Liquid. In PBS, pH 7.2, containing 10mM sodium phosphate, 50mM sodium chloride, and 0.1mM PMSF.

Host Mouse

Immunogen Recombinant human Caspase-7 protein (full length).

Isotype IgG1

Purity Detail Protein G 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.

Source Purified from mouse ascites.

Species Reactivity Hamster, Human, Monkey, Mouse, Rat

UniProt ID P55210

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)