

# 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 »](#)

## Ordering Information

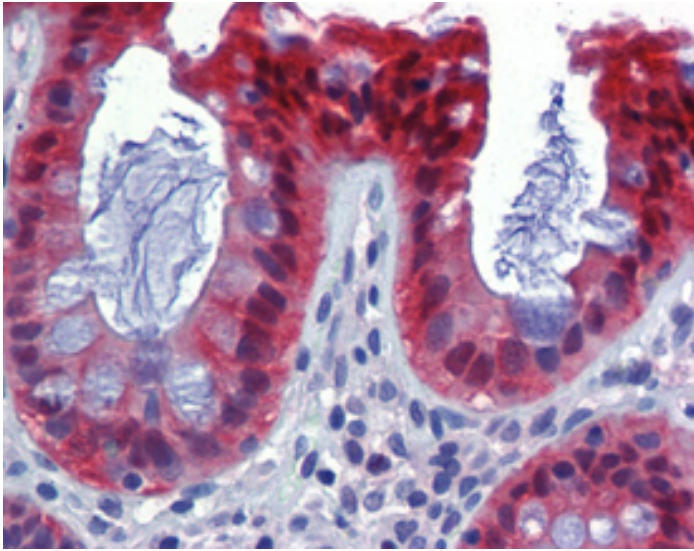
[Order Online »](#)

ADI-AAM-137-C	25µg
ADI-AAM-137-E	100µg

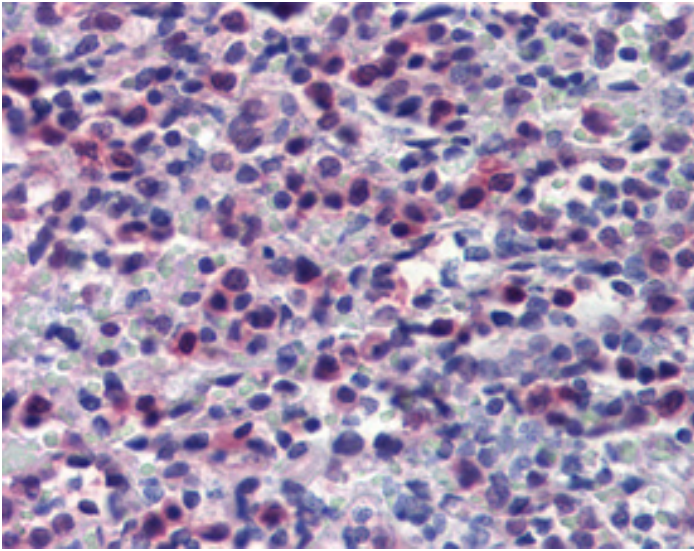
## Manuals, SDS & CofA

[View Online »](#)

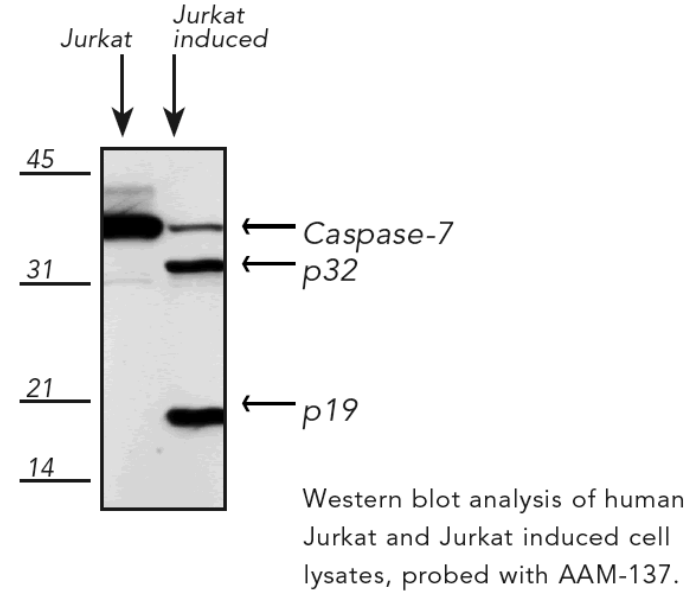




IHC Analysis: Human colon tissue stained with Caspase-7, mAb (10-1-62) at 10µg/ml.



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





<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## 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 <a href="#">Worry-Free Guarantee</a>





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)