

Caspase-3 monoclonal antibody (31A1067)

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

Citations: 43

[View Online »](#)

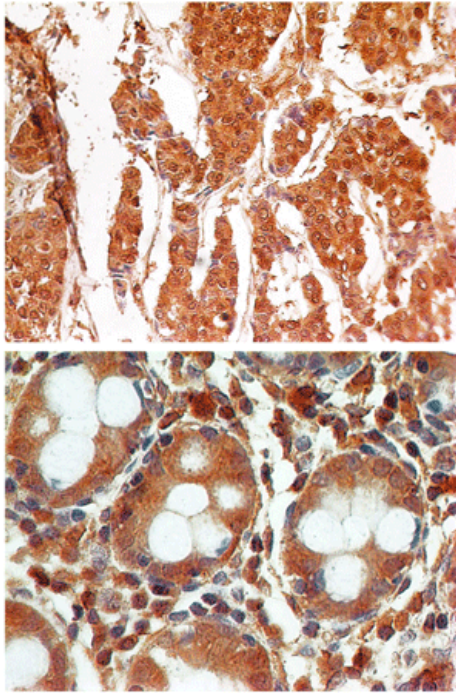
Ordering Information

[Order Online »](#)

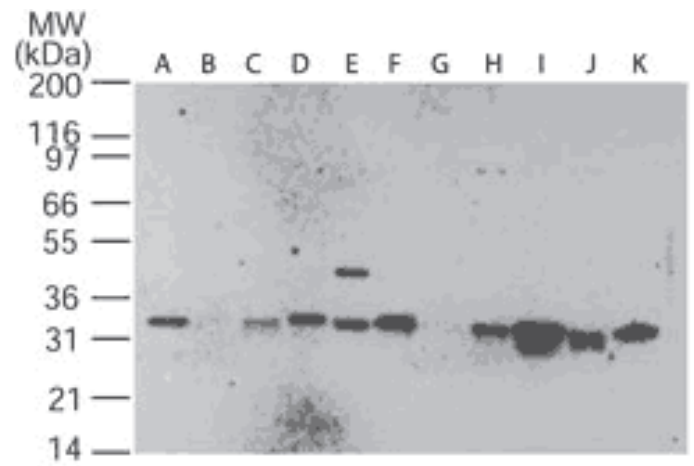
ALX-804-305-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis: Formalin-fixed, paraffin-embedded human breast cancer (top) and normal colon (bottom) stained with Caspase-3, mAb (31A1067) at 4 μ g/ml. Localization can be cytoplasmic and nuclear. Staining in the nucleus is considered to be an indication of active Caspase-3. In most cell types and model systems, cells with active Caspase-3 are undergoing apoptosis.



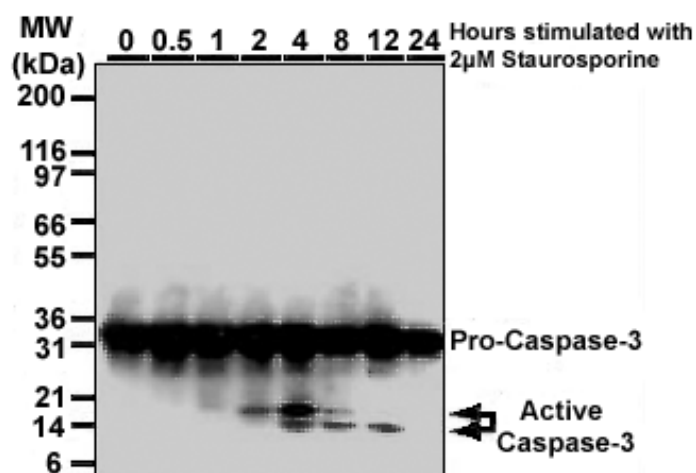


Figure 1: Western blot analysis for detection of Caspase-3 activation in HeLa cells. Cells were treated with 2mM staurosporine for different time periods. Caspase-3 activation is determined by cleavage of procaspase-3, which generates 17 and 12kDa, larger and smaller catalytic subunit, respectively.

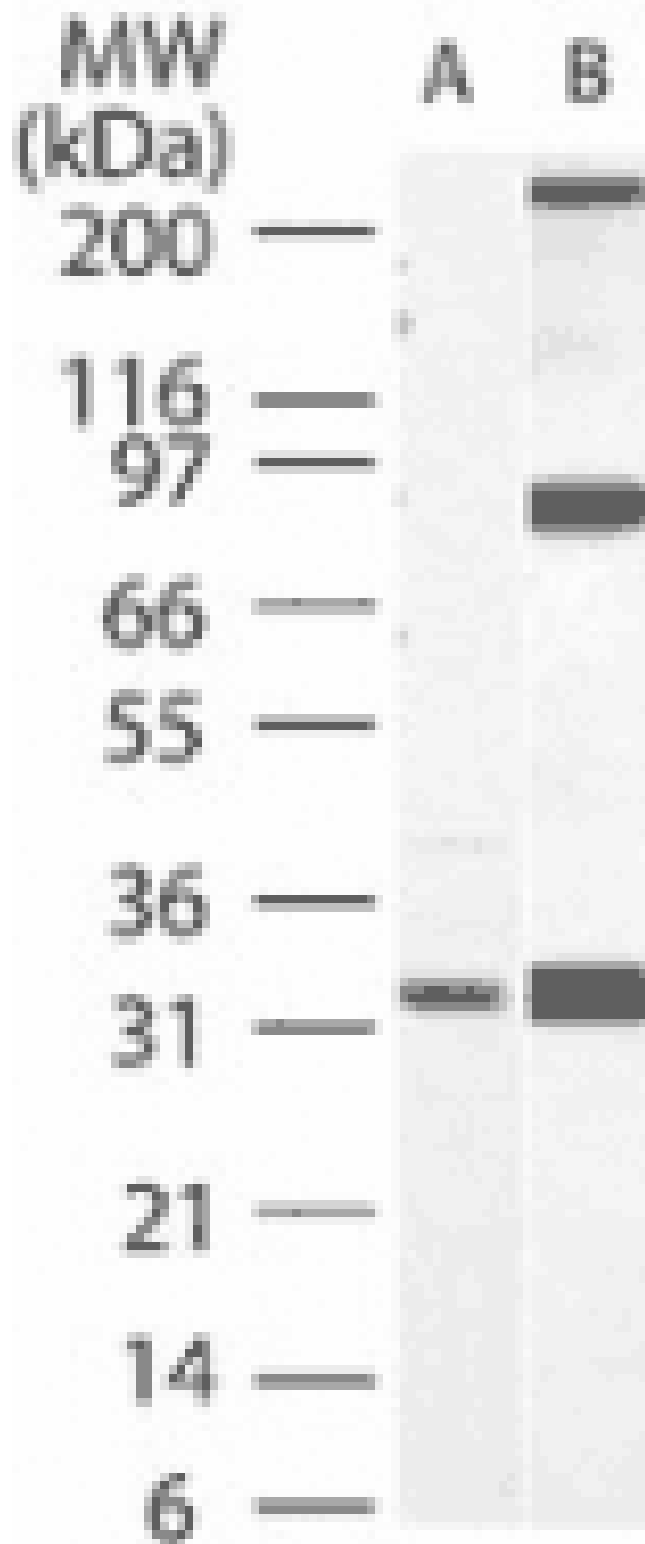


Figure 3: Western blot analysis of Caspase-3 in A) RAW and B) NIH 3T3 using MAb to Caspase-3 (31A1067) (Prod. No. ALX-804-305) 2μg/ml.

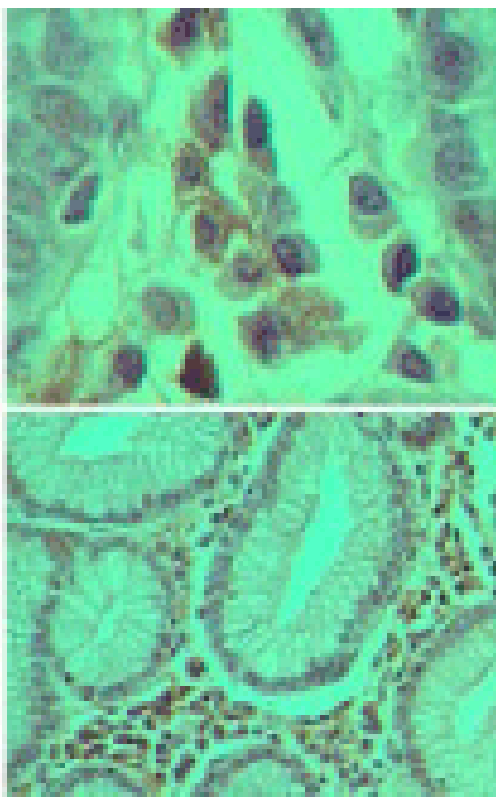


Figure 4: Immunohistochemical analysis of caspase-3 in formalin-fixed, paraffin-embedded human stomach tissue using MAb to Caspase-3 (31A1067) (Prod. No. ALX-804-305) at 5 ug/ml. High (top panel) and low magnification (bottom panel). Caspase-3 staining is cytoplasmic in some cells and nuclear in others. Caspase-3 in the nucleus is considered to be an indication of active caspase 3. In most cell types and model systems, cells with active caspase-3 are undergoing apoptosis.

Handling & Storage

Use/Stability	Stable for 6 months when stored at +4°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	IHC (FS), IHC (PS), WB
Clone	31A1067
Formulation	Liquid. In PBS with 0.05% sodium azide.
Host	Mouse
Immunogen	Recombinant human caspase-3.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Western Blot (2-5µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
Specificity	Recognizes procaspase-3 and cleaved 17kDa catalytic subunit of caspase-3. Recognizes an epitope in the large (17kDa) domain subunit of caspase-3.
UniProt ID	P42574



ENZO LIFE SCIENCES,
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 1 402 44 555
info-fr@enzolifesciences.com

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

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

This antibody is covered by our Worry Free Guarantee