

# BAP31 monoclonal antibody (A1/182)

BAP31 (B cell receptor-associated protein 31) is an integral protein of the ER membrane that forms a large hetero-oligomeric complex with the related BAP29. The cytosolic domain of BAP31 contains two identical caspase recognition sites that are preferentially cleaved by initiator caspases, including caspase-8. After activation of cell surface death receptors, human BAP31 is cleaved into a membrane-embedded fragment called p20, which induces apoptosis when expressed ectopically. The p20 fragment can mediate  $\text{Ca}^{2+}$ -dependent apoptotic cross-talk between the ER and mitochondria, stimulating cytochrome c release. In addition to its role in apoptosis, BAP31 is a component of the ER quality control compartment and functions as a cargo receptor for ER export of transmembrane proteins such as cellubrevin, MHC class I molecules, and cytochrome P450 2C2.

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

Citations: 18

[View Online »](#)

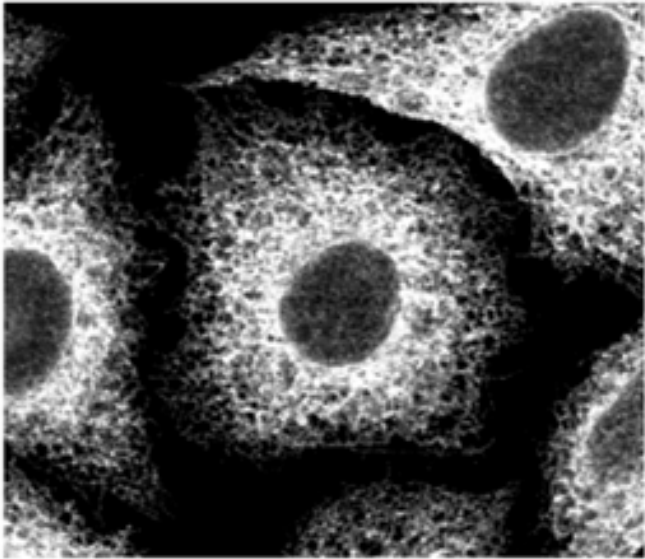
## Ordering Information

[Order Online »](#)

ALX-804-601-C100	100µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



# Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** p28 BAP31, B cell receptor-associated protein BAP31, BCR-associated protein BAP31, CDM, DXS1357E, 6C6-AG tumor-associated antigen, BCAP31

**Application** ELISA, Flow Cytometry, ICC, IHC (FS), IP, WB

**Clone** A1/182

**Formulation** Lyophilized. Protein G-affinity purified and gel filtrated (FPLC).

**Host** Mouse

**Immunogen** BAP31 (BCR associated protein BAP031) from Vero monkey cells.

**Isotype** IgG2a

**Recommendation** Immunocytochemistry (1:1,000)Immunoprecipitation (2µl)Western Blot  
**Dilutions/Conditions** (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Reconstitution** Reconstitute with 100µl PBS containing 0.1% sodium azide.

**Species Reactivity** Human, Monkey

**UniProt ID** P51572 (human)

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



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-  
fr@enzolifesciences.com](mailto:info-fr@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)