

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 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: 19

[View Online »](#)

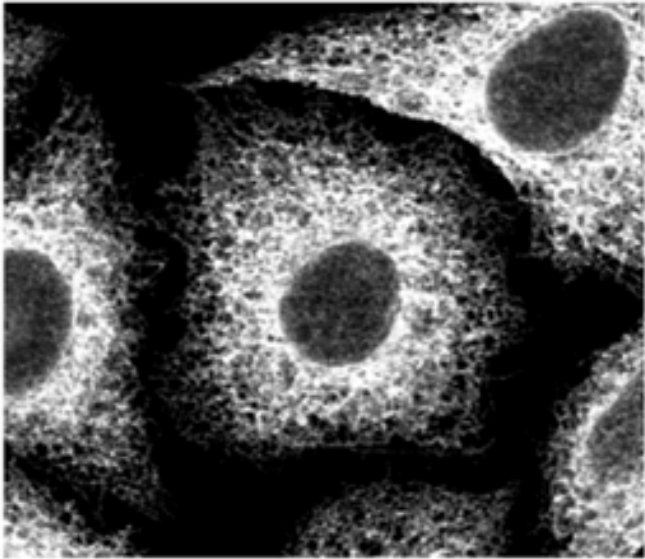
Ordering Information

[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 Dilutions/Conditions	Immunocytochemistry (1:1,000)Immunoprecipitation (2µl)Western Blot (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-
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)