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 Ca²⁺-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

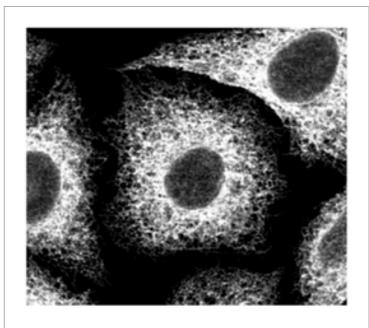
Order Online »

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.

lgG2a

Recommendation Dilutions/Conditions Immunocytochemistry (1:1,000)Immunoprecipitation

(2µI)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

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com