

Bax monoclonal antibody (4F11)

Bax is an approximately 22 kDa pro-apoptotic protein belonging to the Bcl-2 family of proteins, which are characterized by the presence of at least one Bcl-2 homology (BH) domains. Bax normally resides in the cytoplasm, but following an apoptotic stimuli, Bax undergoes a conformational change and translocates to mitochondrial membranes where it inserts and mediates the release of cytochrome c. This activates caspase-3 and perpetuates the apoptotic cascade.

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

Citations: 4

[View Online »](#)

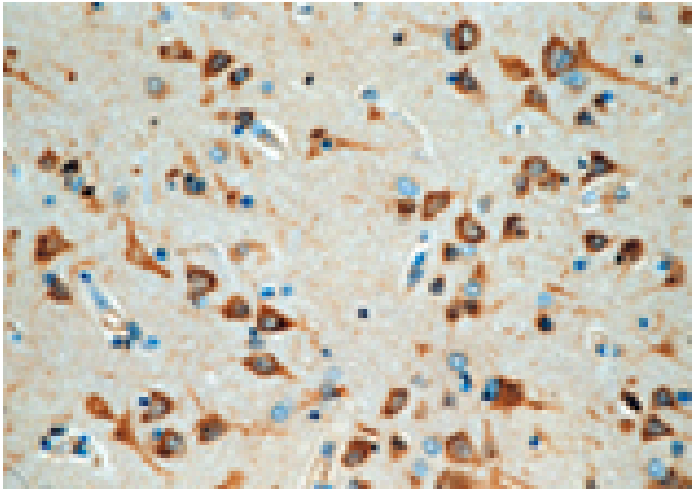
Ordering Information

[Order Online »](#)

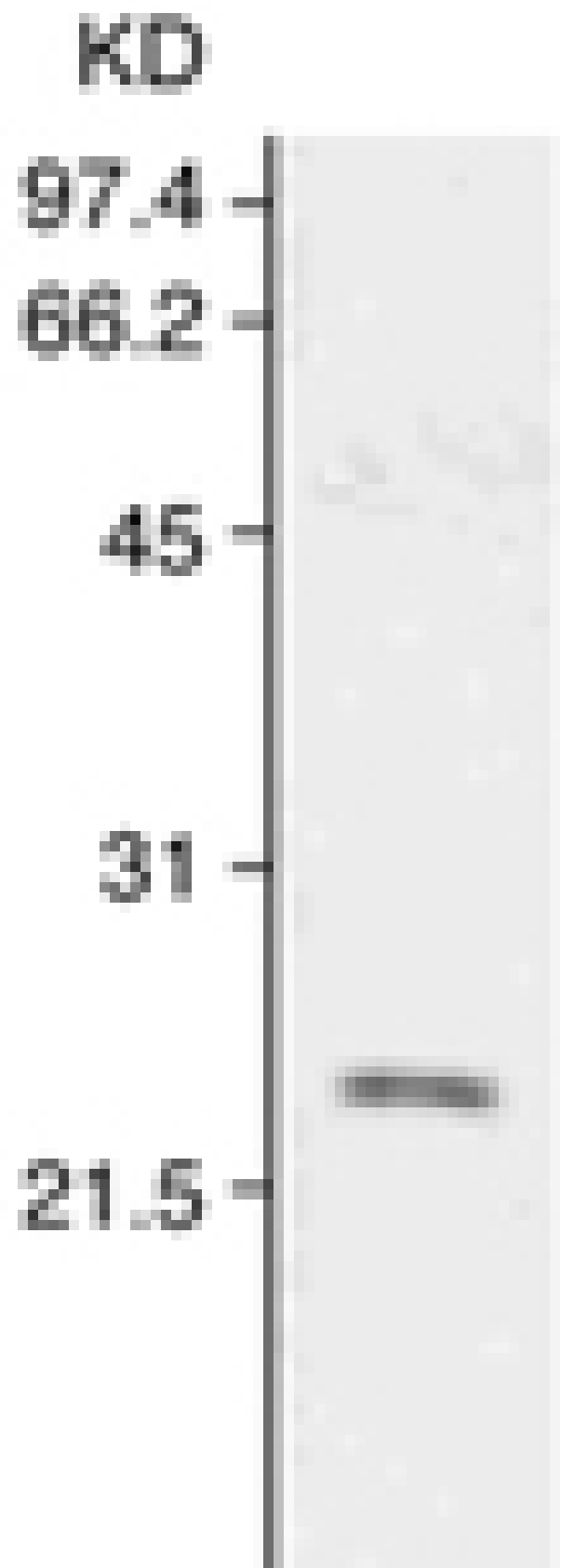
ADI-AAM-140-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of paraffin-embedded section of Alzheimer's diseased (AD) brain, stained with Bax mAb (4F11).



Western blot analysis of human Raji cell lysate, probed with Bax mAb (4F11)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name BCL2L4, BCL2-associated X protein

Application IHC, IP, WB

Application Notes Detects a band of ~22kDa by Western blot.

Clone 4F11

Formulation Lyophilized. In PBS, pH 7.2, containing 1% sucrose. Reconstitute with 100µL of distilled water.

Host Mouse

Immunogen Recombinant human Bax.

Isotype IgG2b

Purity Detail Protein A affinity purified.

Recommendation Immunohistochemistry (10µg/ml) Immunoprecipitation (5-10µg/ml) Western Blot (0.1-1.0µg/ml colorimetric, ECL) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Rat

UniProt ID Q07812

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)