

# 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 »](#)

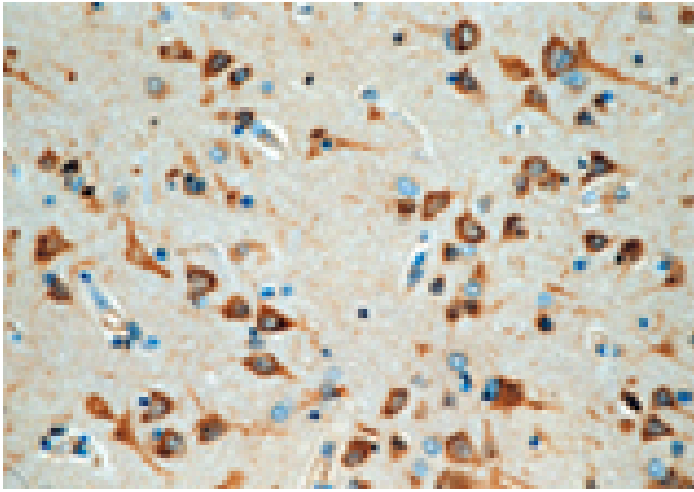
## Ordering Information

[Order Online »](#)

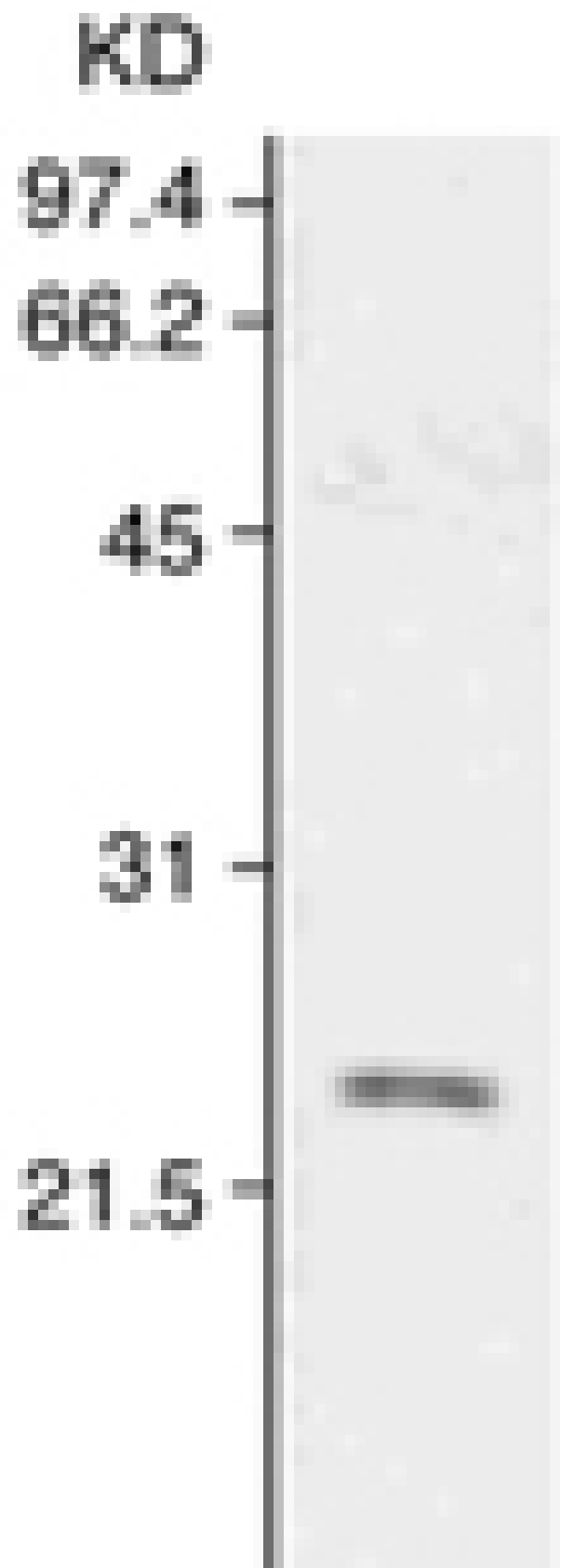
ADI-AAM-140-E	100µg
---------------	-------

## 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 Dilutions/Conditions	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 <a href="#">Worry-Free Guarantee</a> .



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)