

# Bax polyclonal antibody

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: 3 [View Online »](#)

Ordering Information [Order Online »](#)

ADI-AAS-040-D	50µg
ADI-AAS-040-F	200µg

Manuals, SDS & CofA [View Online »](#)

**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)