

# Bak 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](#).

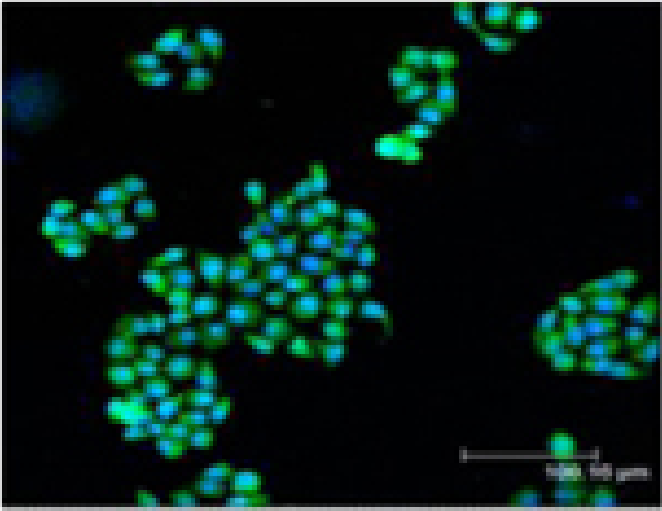
## Ordering Information

[Order Online »](#)

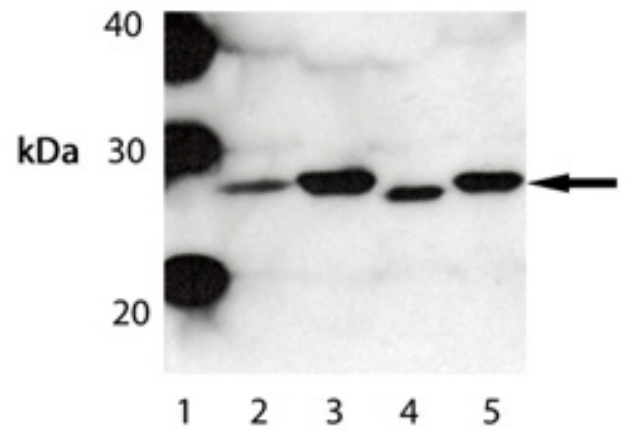
ADI-AAP-030-D	50µg
ADI-AAP-030-F	200µg

## Manuals, SDS & CofA

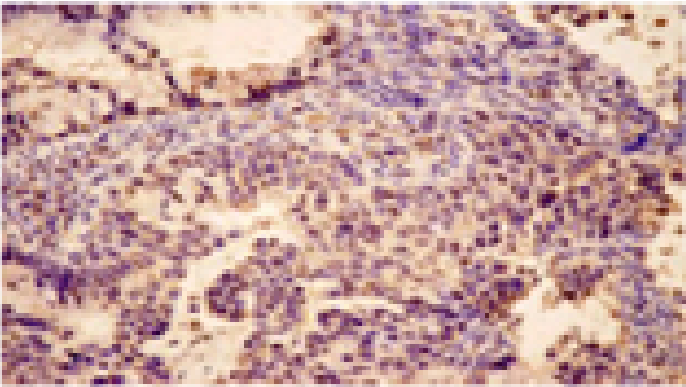
[View Online »](#)



Immunofluorescent analysis (confocal) staining of He La cells using Bak pAb (green); nuclei are stained in blue pseudocolor using DRAQ5.



Western blot analysis of Bak (Prod. no. ADI-AAP-030). Lane 1: MW Marker; Lane 2: A431; Lane 3: CHO-K1; Lane 4: ESK-4; Lane 5: Vero cells.



Immunohistochemistry analysis of human cervical cancer tissue with Bak pAb.

<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Product Details

Alternative Name	BCL2L7
Application	IF, IHC, IP, WB
Application Notes	Detects a band of ~30kDa by Western blot.
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
Gene/Protein Identifier	S58873 (PIR)
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of human Bak.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:50)Immunoprecipitation(1:100)Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Hamster, Human, Monkey, Porcine
UniProt ID	Q16611
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)