

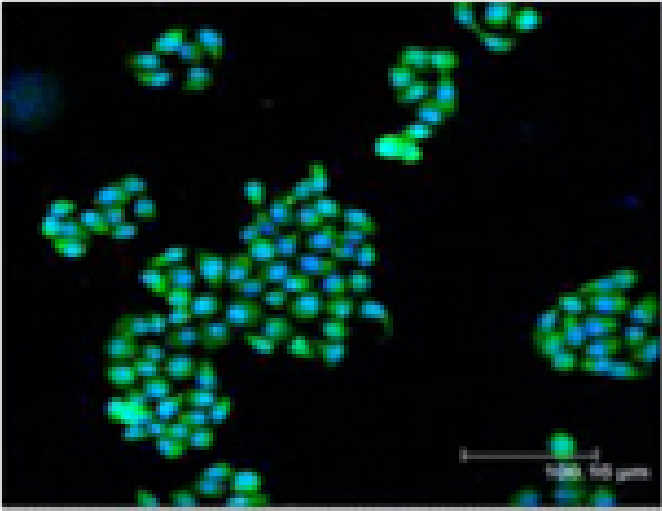
Bax polyclonal antibody

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.

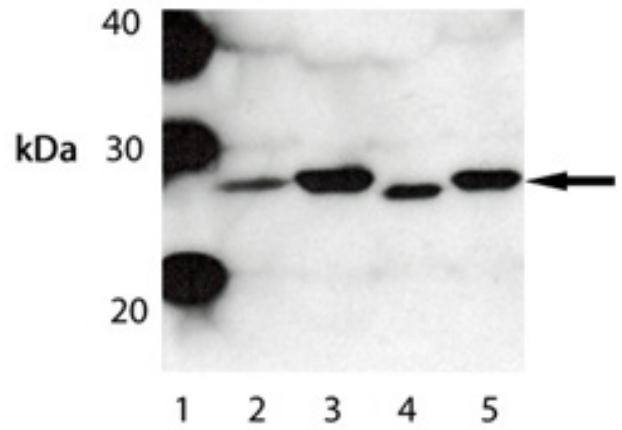
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CoFA

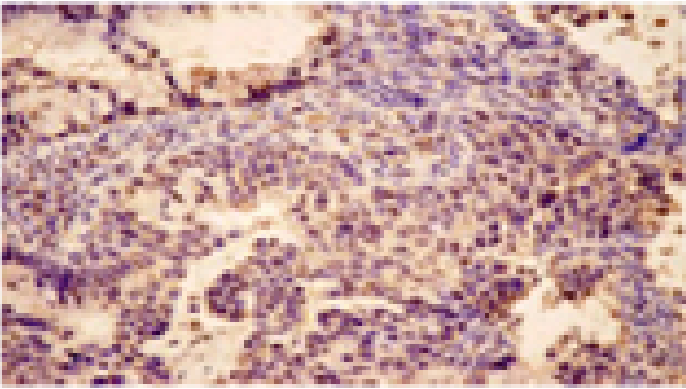
[View Online »](#)



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.

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 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 [Worry-Free Guarantee](#)

Last modified: May 29, 2024



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)