

Bax (active monomer) Recombinant monoclonal antibody (6A7)

Bax and Bak operate in both the endoplasmatic reticulum (ER) and mitochondria as an essential gateway for selected apoptotic signals. A role for Bcl-2 at the ER has been suggested earlier as Bcl-2 may regulate calcium fluxes between the ER and the mitochondria. For a review on Ca^{2+} and its role in apoptosis.

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

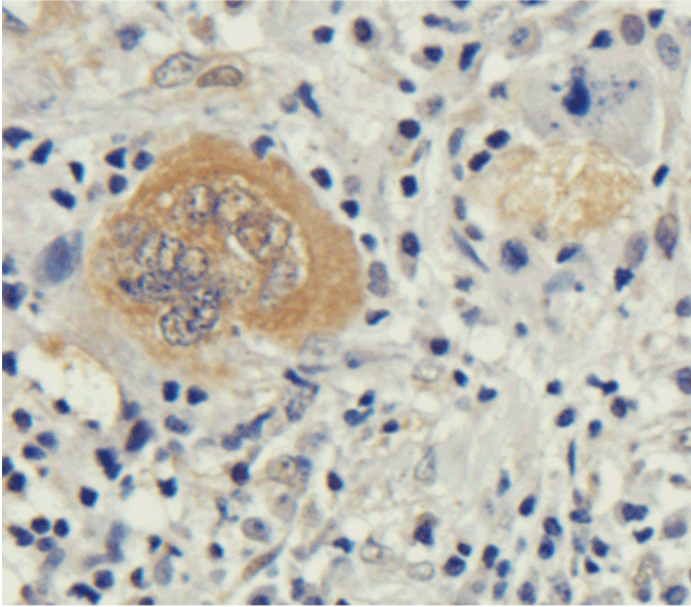
Ordering Information

[Order Online »](#)

ENZ-ABS754-0200	200µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



IHC staining of human lung cancer using anti-Bax antibody. Formalin fixed human lung cancer slices were stained with the antibody at 3 $\mu\text{g/ml}$.

Handling & Storage

Handling	After opening aliquot contents and freeze at -20°C. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	BCL2L4, BCL2-associated X protein
Application	IHC, IP, WB
Clone	6A7
Crossreactivity	In the presence of nonionic detergent, the antibody binds the monomeric form of Bax, but not Bax complexed with either Bcl-XL or Bcl-2. Only recognizes conformationally changed active Bax.
Formulation	Liquid. In PBS with 0.02% Proclin 300.
Host	Mouse
Immunogen	KLH conjugated to peptides corresponding to amino acids 3-16 of murine Bax (CGSGEQLGSGGPTSS), amino acids 3-16 of human Bax (CGSGEQPRGGGPTSS), and amino acids 12-24 shared by Bax from both species (CGPTSSEQIMKTGA).
Isotype	IgG1κ
Purity Detail	Protein A affinity purified.
Species Reactivity	Human, Mouse, Rat
Specificity	Recognizes the active monomeric form of Bax.
UniProt ID	Q07812

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com