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 Ca²⁺ and its role in apoptosis.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

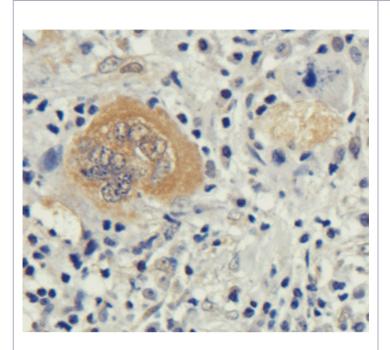
Order Online »

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 were stained with the antibody at 3 μ 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 IgG1ĸ

Purity Detail Protein A affinity purified.

Species Reactivity Human, Mouse, Rat

Specificity Recognizes the active monomeric form of Bax.

UniProt ID Q07812

Last modified: May 29, 2024



info-

eu@enzolifesciences.com