

Bax (NT) antibody (pAb)

Rabbit Anti-Human Bax (Bcl-2 Associated X Protein)

Instruction Manual

Catalog Number	PK-AB718-3351
Synonyms	Bax Antibody:
Description	Apoptosis plays a major role in normal organism development, tissue homeostasis, and removal of damaged cells. Disruption of this process has been implicated in a variety of diseases such as cancer. The Bcl-2 family of proteins is comprised of critical regulators of apoptosis that can be divided into two classes: those that inhibit apoptosis and those that promote cell death. Bax, a pro-apoptotic Bcl-2 family member, is a cytosolic protein that changes conformation and translocates to the mitochondria following apoptotic stimuli. It is thought to share significant functional homology with Bak, another pro-apoptotic Bcl-2 family member, as disruption of bax or bak has little effect on cell death, but mice lacking both genes display multiple developmental defects and cells lacking both show decreased apoptotic capability.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Bax antibody was raised against a peptide corresponding to 16 amino acids near the amino-terminus of human Bax (GenBank accession no. AAA03619).
Purification Method	Ion exchange chromatography purified.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, ICC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually.
Images	Bax antibody can be used for detection of Bax by Western blot at 1 to 4 µg/mL. Antibody can also be used for immunocytochemistry starting at 2 µg/mL. For immunofluorescence start at 2 µg/mL.
References	Lockshin RA, Osborne B, and Zakeri Z. Cell death in the third millennium. Cell Death Differ. 2000; 7:2-7. Cory S, Huang DCS, and Adams JM. The Bcl-2 family: roles in cell survival and oncogenesis. Oncogene 2003; 22:8590-607. Heiser D, Labi V, Erlacher M, et al. The Bcl-2 protein family and its role in the development of neoplastic disease. Exp. Gerontol. 2004; 39:1125-35. Oltvai ZN, Millman CL, and Korsmeyer SJ. Bcl-2 heterodimerizes in vivo with a conserved homolog, Bax, that accelerates programmed cell death. Cell 1993; 74:609-19.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3351P HL-60 Cell Lysate, Cat. No. PK-AB718-1209 Bcl-2 Antibody (NT), Cat. No. PK-AB718-3335 Bak Antibody (NT), Cat. No. PK-AB718-3347

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.