

Bax antibody (pAb)

Rabbit Anti-Human/Mouse/Rat/Monkey Bax (Bcl-2 Associated X Protein)

Instruction Manual

Catalog Number	PK-AB577-3032
Description	The Bcl-2 (B-cell leukemia/lymphoma 2) family proteins function to either inhibit (Bcl-2 and Bcl-XL) or promote (Bax and Bak) apoptosis. Bax contains the conserved Bcl-2 homology 1 (BH1) and 2 (BH2) domains that allow for its homodimerization or heterodimerization with Bcl-2. When Bcl-2 is in excess, apoptosis is inhibited. However, if Bax levels increase in response to a death signal, the cell is pushed toward death.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide corresponding to amino acids near the N-terminus of human Bax.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat, monkey
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit anti-Bax polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 0.01% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	Western blot analysis (0.5-4 µg/ml), immunoprecipitation (10-20 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody detects a 20 kDa protein, corresponding to the apparent molecular weight of Bax on SDS-PAGE immunoblots.

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