

Bcl-xL antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Bcl-xL

Instruction Manual

Catalog Number	PK-AB577-3312
Description	Bcl-xL is a 30 kDa anti-apoptotic protein belonging to the Bcl-2 family. Bcl-xL prevents apoptosis through two different mechanisms: by heterodimerization with an apoptotic protein to inhibit its apoptotic effect, and by its direct pore-forming effect on the outer-membrane of mitochondria to help maintain a normal membrane state under stress conditions. Bcl-xL can be cleaved by caspases upon apoptotic activation, which converts Bcl-xL from an antiapoptotic protein to a proapoptotic cleavage product. The cleavage product of Bcl-xL forms cytochrome c-releasing pores in the mitochondrial membrane to facilitate the apoptosis signaling pathway.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 56 of human Bcl-xL.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) affinity purified, rabbit anti-Bcl-xL 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 blotting (0.5-4 µg/ml) and immunoprecipitation (20 µ/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody detects the 30 kDa Bcl-xL protein. It does not react with other Bcl-2 family members.

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