

Instruction Manual

Catalog Number	PK-AB577-3172
Description	Bid, a BH3 domain-containing proapoptotic Bcl-2 family member, is localized in the cytosolic fraction of cells as an inactive precursor. Its active form is generated upon proteolytic cleavage by caspase-8 in the Fas signaling pathway. Cleaved Bid translocates to mitochondria and releases its potent proapoptotic activity, which in turn induces cytochrome c release and mitochondrial damage. The cytochrome c releasing activity of Bid was antagonized by Bcl-2. Mutation in the SH3 domain can diminish the cytochrome c releasing activity. In the animal model studies, Bid-deficient mice are found resistant to the lethal effects of death factor signals relayed through Fas.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 37 of human Bid.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit anti-Bid 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 (5-10 µg/ml), and immunohistochemistry (20-40 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody detects 22 kDa human Bid.

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