

Bid (IN) antibody (pAb)

Rabbit Anti-Human/Mouse Bid (Member of the Bcl-2 Family)

Instruction Manual

Catalog Number	PK-AB718-3355
Synonyms	Bid 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. Bid, a pro-apoptotic Bcl-2 family member, is cleaved by caspase-8 in response to apoptotic signals, exposing the Bcl-2 homology 3 (BH3) domain which is normally buried in the full-length protein. The cleaved complex is myristoylated and translocated to the mitochondrial membrane where it may induce mitochondrial Bax and Bak to oligomerize.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Bid antibody was raised against a peptide corresponding to 14 amino acids near the middle of human Bid (GenBank accession no. AAH36364).
Purification Method	Ion exchange chromatography purified.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
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 INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Application have to be determined individually. Bid antibody can be used for detection of Bid by Western blot at 0.5 to 2 µg/mL.
Images	Available upon request.
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. Wang K, Yin XM, Chao DT, et al. BID: a novel BH3 domain-only death agonist. Genes Dev. 1996; 10:2859-69.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3355P A549 Cell Lysate, Cat. No. PK-AB718-1203 Bcl-2 Antibody (NT), Cat. No. PK-AB718-3335 Bak Antibody (NT), Cat. No. PK-AB718-3347 Bax Antibody (NT), Cat. No. PK-AB718-3351 Bid Antibody (CT), Cat. No. PK-AB718-3353

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