

Bim / BOD (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Bim (Bcl-2 Interacting Mediator of Cell Death)

Instruction Manual

Catalog Number	PK-AB718-2065
Synonyms	Bim Antibody: BOD
Description	Members in the Bcl-2 family are critical regulators of apoptosis by either inhibiting or promoting cell death. Bcl-2 homology 3 (BH3) domain is a potent death domain. BH3 domain containing proapoptotic proteins, including Bad, Bax, Bid, Bik, and Hrk, form a growing subclass of the Bcl-2 family. A novel BH3 domain containing protein was recently identified and designated Bim or BOD in human, mouse and rat. Bim/BOD interacts with diverse members in the pro-survival Bcl-2 subfamily including Bcl-2, Bcl-xL and Bcl-w. Bim/BOD induces apoptosis. The messenger RNA of Bim is ubiquitously expressed in multiple tissues and cell lines.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Bim antibody was raised against a peptide corresponding to amino acids near the center of human Bim. The sequence is identical to that of mouse and differs from that of rat by one amino acid.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
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, IHC Note: 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. Bim antibody can be used for detection of Bim by Western blot at 1 µg/mL. A 23 kDa band can be detected. Antibody can also be used for immunohistochemistry starting at 20 µg/mL.
Images	Available upon request.
References	O'Connor L, Strasser A, O'Reilly LA, et al. Bim: a novel member of the Bcl-2 family that promotes apoptosis. EMBO J. 1998; 17:384-395. Hsu SY, Lin P, and Hsueh AJ BOD (Bcl-2-related ovarian death gene) is an ovarian BH3 domain-containing proapoptotic Bcl-2 protein capable of dimerization with diverse antiapoptotic Bcl-2 members. Mol. Endocrinol. 1998; 12:1432-40.
Images	Available upon request.
Related Products	K562 Cell Lysate, Cat. No. PK-AB718-1204 Bim Blocking Peptide, Cat. No. PK-AB718-2065P

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