

Bfl-1 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Bfl-1 (Bcl-2 Related Protein A1)

Instruction Manual

Catalog Number	PK-AB718-3875
Synonyms	Bfl-1 Antibody: Bcl-2-related protein A1, Hemopoietic-specific early response protein
Description	Apoptosis plays a major role in normal organism development, tissue homeostasis, and removal of damaged cells and is caused by caspase activation. Proteins that comprise the Bcl-2 family appear to control the activation of these enzymes. One such member is multi-domain antiapoptotic protein Bfl-1, which is overexpressed in stomach and other cancers. Bfl-1 can interact with Bax and suppress apoptosis by inhibiting the release of cytochrome c and caspase-3 activation. It is upregulated in cisplatin-resistant human bladder tumors, suggesting that its expression may be important for cisplatin resistance and inhibition of apoptosis in cancer cells. At least two isoforms of Bfl-1 are known to exist. Presumably due to post-translational modifications, Bfl-1 is often observed at higher molecular weight in SDS-PAGE than its predicted molecular weight would suggest.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Bfl-1 antibody was raised against a 14 amino acid peptide from near the amino terminus of human Bfl-1.
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, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Bfl-1 antibody can be used for the detection of Bfl-1 by Western blot at 1 - 2 µg/mL. Presumably due to post-translational modifications, Bfl-1 is often observed at higher molecular weight in SDS-PAGE than its predicted molecular weight would suggest. For immunofluorescence start at 20 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. Choi SS, Park IC, Yun JW, et al. A novel Bcl-2 related gene, Bfl-1, is overexpressed in stomach cancer and preferentially expressed in bone marrow. Oncogene 1995; 11:1693-8. Kim JK, Kim KD, Lee E, et al. Up-regulation of Bfl-1/A1 via NF-κB activation in cisplatin-resistant human bladder cancer cell line. Cancer Lett. 2004; 212:61-70. Zhang H, Cowan-Jacob SW, Simonen M, et al. Structural basis of BFL-1 for its interaction with BAX and its anti-apoptotic action in mammalian and yeast cells. J. Biol. Chem. 2000; 275:11092-9.
Images	Available upon request.
Related Products	Kidney Lysate, Cat. No. PK-AB718-1405 Bcl-2 Antibody, Cat. No. PK-AB718-3335; Bax Antibody, Cat. No. PK-AB718-3351 Bfl-1 Antibody, Cat. No. PK-AB718-3873
FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.	