

# Bnip3L / Nix (IN) antibody (pAb)

Rabbit Anti-Human Bnip3L (Member of the Bcl-2 Family)

## Instruction Manual

Catalog Number	PK-AB718-2289
Synonyms	Bnip3L Antibody: Nix
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 pro-apoptotic proteins, including Bad, Bax, Bid, Bik, Hrk, Nip3, and Bim, form a growing subclass of the Bcl-2 family. A novel BH3 domain containing protein was recently identified and designated Bnip3L, Bnip3 $\alpha$ , and Nix (for Nip3-like protein X). Bnip3L/Bnip3 $\alpha$ /Nix is a homolog of the E1B 19K/Bcl-2 binding and pro-apoptotic protein Bnip3. Overexpression of Bnip3L induces apoptosis. Bnip3L interacts with and overcomes suppresses by Bcl-2 and Bcl-xL. Bnip3L is localized in mitochondria. The messenger RNA of Bnip3L is ubiquitously expressed in human tissues. Bnip3L and Bnip3 form a new subfamily of the pro-apoptotic mitochondrial proteins.
Quantity	100 $\mu$ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Bnip3L antibody was raised against a peptide (DAQHESGQSSSRGSSH) corresponding to amino acids 77 to 92 of human origin, which are identical to those of mouse Bnip3L .
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
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 $\mu$ l 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 <span style="float:right">INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually.</span> Bnip3L antibody can be used for detection of Bnip3L by Western blot at 1 $\mu$ g/mL. The approximate 40 kDa top band represents monomer of Bnip3L (4). Antibody can also be used for immunohistochemistry starting at 2 $\mu$ g/mL. For immunofluorescence start at 10 $\mu$ g/mL.
Images	Available upon request.
References	Matsushima M, Fujiwara T, Takahashi E, et al. Isolation, mapping, and functional analysis of a novel human cDNA (BNIP3L) encoding a protein homologous to human NIP3. <i>Genes Chromosomes Cancer</i> 1998;21:230-5 Yasuda M, Han JW, Dionne CA, Boyd JM, Chinnadurai G. BNIP3 $\alpha$ : a human homolog of mitochondrial proapoptotic protein BNIP3. <i>Cancer Res</i> 1999;59:533-7 Chen G, Cizeau J, Vande Velde C, et al. Nix and Nip3 form a subfamily of pro-apoptotic mitochondrial proteins. <i>J Biol Chem</i> 1999;274(1):7-10 Imazu T, Shimizu S, Tagami S, et al. Bcl-2/E1B 19 kDa-interacting protein 3-like protein (Bnip3L) interacts with bcl-2/Bcl-xL and induces apoptosis by altering mitochondrial membrane permeability. <i>Oncogene</i> . 1999;18:4523-9 (RD1200)
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-2289P K562 Cell Lysate, Cat. No. PK-AB718-1204

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