

Bcl-2 (IN) antibody (pAb)

Rabbit Anti-Human B-cell Lymphoma/Leukemia 2

Instruction Manual

Catalog Number	PK-AB718-3337
Synonyms	Bcl-2 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. Bcl-2 is the founding member of a family of over 20 proteins that are critical regulators of apoptosis. These can be divided into two classes: those that inhibit apoptosis and those that promote cell death. Bcl-2 is an inner mitochondrial membrane protein that inhibits apoptosis. It is thought to act by interacting with pro-apoptotic Bcl-2 family members such as Bak and Bad. Overexpression of Bcl-2 has been linked to human cancers such as B-cell lymphoma and prostate cancer.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Bcl-2 antibody was raised against a peptide corresponding to 14 amino acids near the middle of human Bcl-2 (GenBank accession no. AAH27258).
Purification Method	Ion exchange chromatography purified.
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 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, ICC 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. Bcl-2 antibody can be used for detection of Bcl-2 by Western blot at 1 to 4 µg/mL. Antibody can also be used for immunocytochemistry starting at 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. Hockenbery D, Nunez G, Milliman C, et al. Bcl-2 is an inner mitochondrial membrane protein that blocks programmed cell death. Nature 1990; 348:334-6.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3337P A-20 Cell Lysate, Cat. No. PK-AB718-1288 Bcl-2 Antibody (NT), Cat. No. PK-AB718-3335 Bad Antibody (CT), Cat. No. PK-AB718-3343 Bak Antibody (NT), Cat. No. PK-AB718-3347

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