

# Bak (NT) antibody (pAb)

## Rabbit Anti-Human/Mouse Bak (Bcl-2 Antagonist/Killer)

### Instruction Manual

Catalog Number	PK-AB718-3347
Synonyms	Bak 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. Bak, a pro-apoptotic Bcl-2 family member, is an oligomeric protein that localizes to the mitochondria. It is thought to share significant functional homology with Bax, another pro-apoptotic Bcl-2 family member, as disruption of bak or bax has little effect on cell death, but mice lacking both genes display multiple developmental defects and cells lacking bak and bax show decreased apoptotic capability.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Bak antibody was raised against a peptide corresponding to 13 amino acids near the amino-terminus of human Bak (GenBank accession no. Q16611).
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, 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. Bak antibody can be used for detection of Bak by Western blot at 1 to 2 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. Kiefer M, Brauer MJ, Powers VC, et al. Modulation of apoptosis by the widely distributed Bcl-2 homologue Bak. Nature 1995; 374:736-9.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3347P L1210 Cell Lysate, Cat. No. PK-AB718-1284 Bcl-2 Antibody (NT), Cat. No. PK-AB718-3335 Bcl-2 Antibody (IN), Cat. No. PK-AB718-3337 Bax Antibody (NT), Cat. No. PK-AB718-3351
FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.	