

# Bcl-B / Bcl-2L10 antibody (pAb)

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

### Instruction Manual

Catalog Number	PK-AB718-3163
Synonyms	Bcl-B Antibody: Boo, Diva, Nrh
Description	Members in the Bcl-2 family are critical regulators of apoptosis by either inhibiting or promoting cell death. Bcl-B is a recently discovered anti-apoptotic member of the Bcl-2. Unlike the mouse homolog (also known as Diva/Boo) which is predominantly expressed in ovary and testis, the human Bcl-B protein is widely expressed. Also, the human Bcl-B functions by binding to and suppressing the apoptotic activity of Bax, whereas the mouse homolog binds Bak and also interacts with the apoptosis protein Apaf-1. Despite its predicted molecular weight, Bcl-B is often at higher molecular weights, presumably due to post-translational modifications.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Bcl-B antibody was raised against a 16 amino acid peptide from near the amino terminus of human Bcl-B.
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 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 <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>
Images	Bcl-B antibody can be used for the detection of Bcl-B by Western blot at 1 µg/mL. Despite its predicted molecular weight, Bcl-B is often at higher molecular weights, presumably due to post-translational modifications. Antibody can also be used for immunocytochemistry starting at 10 µg/mL.
References	Cory S, Huang DCS, and Adams JM. The Bcl-2 family: roles in cell survival and oncogenesis. <i>Oncogene</i> 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. <i>Exp. Gerontol.</i> 2004; 39:1125-35. Ke N, Godzik A, and Reed JC. Bcl-B: A novel Bcl-2 family member that differentially binds and regulates Bax and Bak. <i>J. Biol. Chem.</i> 2001; 276:12481-4. Song Q, Kuang Y, Dixit VM, et al. Boo, a negative regulator of cell death, interacts with Apaf-1. <i>EMBO J.</i> 1999; 18:167-78.
Images	Available upon request.
Related Products	Apaf-1 Antibody, Cat. No. PK-AB718-2015 Bak Antibody, Cat. No. PK-AB718-3347 Bax Antibody, Cat. No. PK-AB718-3351 Bcl-2Antibody, Cat. No. PK-AB718-3335 Bcl-B Peptide, Cat. No. PK-AB718-3163P Jurkat Lysate, Cat. No. PK-AB718-1205

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