

Rabbit Anti-Human/Mouse/Rat B-cell Lymphoma/Leukemia 10 (Involved in Apoptosis and Tumor Genesis)

Instruction Manual

Catalog Number	PK-AB718-2161
Synonyms	Bcl-10 Antibody: CIPER, CLAP
Description	Apoptosis is related to many diseases including cancer. Cell death signals are transduced by death domain (DD) and caspase recruitment domain (CARD) containing molecules and a caspase family of proteases. CARD containing cell death regulators include ARC, RAIDD, Apaf-1, caspase-9, and caspase-2. A novel CARD containing protein was recently identified by several groups and designated Bcl10, CIPER, mE10, CARMEN, CLAP. Bcl10 is a cellular homolog of the equine herpesvirus-2 E-10 gene. Overexpression of Bcl10 induces JNK, p38, and NF-κB activation. Bcl10 interacts with caspase-9 and enhances pro-caspase-9 processing and induces apoptosis through caspase-9 activation. Bcl10 exhibits a variety of mutations in MALT lymphomas and in B and T cell lineage lymphomas indicating that it may be commonly involved in the pathogenesis of human malignancy. Bcl10 is expressed in many human and murine tissues and cell lines.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Bcl10 antibody was raised against a peptide corresponding to amino acids near the amino terminus of human Bcl10.
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, ICC, IF 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-10 antibody can be used for detection of BCL10 by Western blot at 0.5 µg/mL. dilution. An approximately 31 kDa band can be detected. Antibody can also be used for immunocytochemistry starting at 1 µg/mL. For immunofluorescence start at 10 µg/mL.
Images	Available upon request.
References	Willis TG, Jadayel DM, Du MQ, et al. Bcl10 is involved in t(1;14)(p22;q32) of MALT B cell lymphoma and mutated in multiple tumor types. <i>Cell</i> 1999;96(1):35-45 Koseki T, Inohara N, Chen S, et al. CIPER, a novel NF κB-activating protein containing a caspase recruitment domain with homology to Herpesvirus-2 protein E10. <i>J Biol Chem</i> 1999;274(15):9955-61 Yan M, Lee J, Schilbach S, Goddard A, Dixit V. mE10, a novel caspase recruitment domain-containing proapoptotic molecule. <i>J Biol Chem</i> 1999;274(15):10287-92 Thome M, Martinon F, Hofmann K, et al. Equine herpesvirus-2 E10 gene product, but not its cellular homologue, activates NF-κB transcription factor and c-Jun N-terminal kinase. <i>J Biol Chem</i> 1999;274(15):9962-8
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-2161P; Raji Cell Lysate, Cat. No. PK-AB718-1207 Apaf-1 Antibody (NT), Cat. No. PK-AB718-2013; ARC Antibody (CT), Cat. No. PK-AB718-2081 Caspase-9 Antibody (IN2), Cat. No. PK-AB718-2073

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