

## Bcl-10

### B-cell lymphoma 10

human, recombinant, Sf9 insect cells

Cat. No.	Amount
PR-769	6 µg

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -80°C

### Avoid freeze / thaw cycles

#### Form

Liquid. Supplied in 20 mM Tris-HCl pH 8.0, 100 mM KCl, 0.2 mM EDTA, 1 mM DTT and 20% glycerol.

#### Activity

100 ng are required for a protein-protein interaction assay.

#### Application

Recombinant Bcl-10 protein can be used for

- 1) protein-protein interaction assay
- 2) for *in vitro* transcription assay
- 3) for cell growth assay.

#### Molecular Weight

33 kDa

#### Purity

> 95% by SDS-PAGE

### Description

Bcl-10 is an intracellular signaling protein identified from the t (p22;q32) breakpoint in MALT lymphomas. The human Bcl-10 gene encodes a protein of 233 amino acids containing an N-terminal Caspase Recruitment Domain (CARD) that mediates self-oligomerization and the C-terminal region of Bcl-10. Bcl-10 is expressed ubiquitously with high expression levels in lymphoid tissues and in the developing central nerve system. Mutations have been found in cases of follicular lymphoma and diffuse large B cell lymphoma. Bcl-10 is a positive regulator of antigen-receptor induced NF-κB activation and Bcl-10-initiated activation of NF-κB can be inhibited by cotransfection of dominant-negative mutants of TRAF2, NIK, IKKα or IκBα. The wild type human Bcl-10 protein was expressed in baculovirus system and purified by an affinity column in combination with FPLC chromatography. Purified protein is greater than 95% homogeneous and contains no detectable proteases, DNase, and RNase activity.

### Selected References:

- Willis *et al.* (1999) Bcl10 is involved in t(1;14)(p22;q32) of MALT B cell lymphoma and mutated in multiple tumor types. *Cell* **96**:35.
- Srinivasula *et al.* (1999) CLAP, a novel caspase recruitment domain-containing protein in the tumor necrosis factor receptor pathway, regulates NF-kappaB activation and apoptosis. *J. Biol. Chem.* **274**:17946.
- Costanzo *et al.* (1999) c-E10 is a caspase-recruiting domain-containing protein that interacts with components of death receptors signaling pathway and activates nuclear factor-kappaB. *J. Biol. Chem.* **274**:20127.
- Koseki *et al.* (1999) CIPER, a novel NF kappaB-activating protein containing a caspase recruitment domain with homology to Herpesvirus-2 protein E10. *J. Biol. Chem.* **274**:9955.
- Thome *et al.* (1999) Equine herpesvirus-2 E10 gene product, but not its cellular homologue, activates NF-kappaB transcription factor and c-Jun N-terminal kinase. *J. Biol. Chem.* **274**:9962.
- Ruland *et al.* (2001) Bcl10 is a positive regulator of antigen receptor-induced activation of NF-kappaB and neural tube closure. *Cell* **104**:33.
- Rebeaud *et al.* (2008) The proteolytic activity of the paracaspase MALT1 is key in T cell activation. *Nature Immunology* **9**:272-281.