

## Bcl10 antibody (mAb)

**Catalog Nos:** 39393, 39394

**RRID:** AB\_2615055

**Clone:** 151

**Isotype:** IgG

**Application(s):** IP, WB

**Reactivity:** Human

**Quantities:** 200 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 33 kDa

**Background:** Bcl10 (B-cell CLL/lymphoma 10) protein promotes apoptosis and activation of NFκB (NFκB p50 & NFκB p65) via NIK and IKK (IKKα, IKKβ, IKKi/IKKε and IKKγ). Bcl10 protein contains a caspase recruitment domain (CARD) which interacts with other CARD domain containing proteins to regulate NFκB signaling. It is also a substrate for MALT1.

**Immunogen:** This Bcl10 antibody was raised against full-length recombinant human Bcl10 protein.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

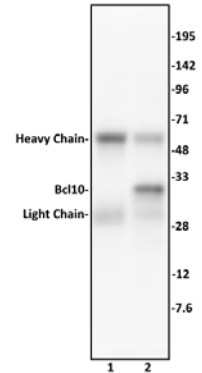
Applications Validated by Active Motif:

IP: 10 µg per IP

WB: 0.5 - 2 µg/ml dilution

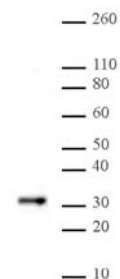
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



### Bcl10 antibody (mAb) (Clone 151) tested by Immunoprecipitation.

10 µg of Bcl10 antibody was used to immunoprecipitate Bcl10 from 500 µg of K562 whole cell extract (lane 2). 10 µg of mouse IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the Bcl10 antibody at a 1 µg/ml dilution.



### Bcl10 antibody (mAb) (Clone 151) tested by Western blot.

Bcl10 detection by Western blot. The analysis was performed using 20 µg HeLa nuclear extract and Bcl10 mAb (Clone 151) at a 1 µg/ml dilution.