

I $\kappa$ B $\alpha$  antibody (mAb)**Catalog No:** 40903**RRID:** AB\_2753150**Clone:** 6A920**Application(s):** WB**Reactivity:** Human, Mouse**Quantity:** 100  $\mu$ g**Purification:** Affinity Purified**Host:** Mouse**Isotype:** IgG1**Concentration:** 1  $\mu$ g/ $\mu$ l**Molecular Weight:** 40 kDa

**Background:** I $\kappa$ B $\alpha$  – NF $\kappa$ B (NF $\kappa$ B p50 & NF $\kappa$ B p65) signaling is controlled to a large extent by the sequestration of the NF $\kappa$ B complex in the cytoplasm by its association with one of the I $\kappa$ B family of proteins. I $\kappa$ B $\alpha$  is phosphorylated at Ser32 and Ser36 (I $\kappa$ B $\alpha$  phospho Ser32,36) by the I $\kappa$ B Kinase (IKK) complex, resulting in the degradation of I $\kappa$ B and the nuclear translocation of NF $\kappa$ B.

**Immunogen:** This I $\kappa$ B $\alpha$  antibody was raised against a recombinant protein corresponding to amino acid residues 32-291 of human I $\kappa$ B $\alpha$ .

**Buffer:** PBS containing 0.02% sodium azide. Sodium azide is highly toxic.

**Application Notes:**

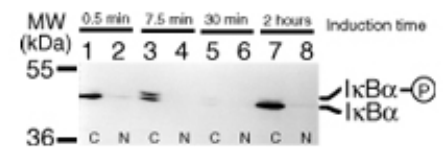
Applications Validated by Active Motif:

WB: 1 - 2  $\mu$ g/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. 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.

**I $\kappa$ B $\alpha$  mAb tested by Western blot.**

The anti-I $\kappa$ B $\alpha$  detects both phosphorylated and non-phosphorylated forms of I $\kappa$ B $\alpha$  by Western blot. Jurkat cells ( $\sim 1 \times 10^7$ ) were treated for indicated time periods with 10 nm/ml PMA, 1  $\mu$ M ionomycin and anti-CD28 (1:10,000 dilution). 10  $\mu$ l of cytoplasmic (C) and nuclear extract (N) were resolved on a 8.75% SDS-PAGE and transferred to Immobilon membrane. Figure is courtesy of Dr. Shao-Cong Sun at Penn State Univ. College of Medicine.