

IκBα phospho Ser32,36 antibody (mAb)

Catalog No: 40904

RRID: AB_2793427

Clone: 39A1431

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Isotype: IgG1

Concentration: 1 µg/µl

Molecular Weight: 40 kDa

Background: IκBα phospho Ser32,36 – NFκB (NFκB p50 & NFκB p65) signaling is controlled to a large extent by the sequestration of the NFκB complex in the cytoplasm by its association with one of the IκB family of proteins. IκBα is phosphorylated at Ser32 and Ser36 by the IκB Kinase (IKK) complex, resulting in the degradation of IκB and the nuclear translocation of NFκB.

Immunogen: This IκBα phospho Ser32,36 antibody was raised against a synthetic peptide containing phosphorylated serines at amino acid residues 32 and 36 of human IκBα.

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

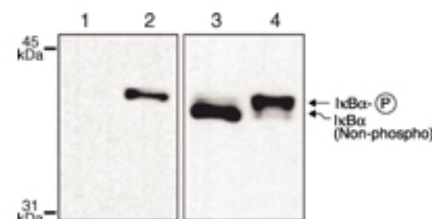
Application Notes:

Applications Validated by Active Motif:

WB: 1 - 2 µg/ml dilution

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.



IκBα phospho Ser32,36 mAb tested by Western blot.

Jurkat cells were treated for 30 minutes with 100 µg/ml ALLN (N-Acetyl-Leu-Leu-Norleucinal), a calpain and proteasome inhibitor that prevents IκBα dephosphorylation, followed by incubation with (lanes 2 & 4) or without 1 nM TNF-α (1 & 3). The membranes were blotted with the anti-phospho-IκBα (lanes 1 & 2) or anti-IκBα, which recognizes both phosphorylated and non-phosphorylated forms of IκBα. The data shows that IκBα phospho Ser32,36 mAb detects specifically the phosphorylated forms of IκBα.