

NFκB p50 antibody (pAb)

Catalog No: 39037

RRID: AB_2793150

Isotype: IgG

Application(s): WB

Reactivity: Human, Mouse

Volume: 200 μl

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.1 μg/μl

Molecular Weight: 50 kDa

Background: NFκB p50 is a member of the NFκB family of transcription factors that are involved in processes such as inflammation, immunity, differentiation, cell growth, tumorigenesis and apoptosis. The Rel family includes c-Rel, RelA (NFκB p65, NFκB3) and RelB (I-Rel). Rel proteins form dimers with the NFκB p50 (proteolytically processed from p105/NFκB1) and p52 (proteolytically processed from p100/NFκB2) members of the NFκB family. The NFκB p65/p50 heterodimer is the most abundant form of NFκB.

NFκB signaling is negatively regulated by the sequestration of the NFκB complex in the cytoplasm by its association with the IκB family of inhibitory proteins. The IκB Kinase (IKK) complex is the key enzyme involved in the activation and translocation of NFκB. The IKK complex is composed of two catalytic subunits (IKKα and IKKβ) and a regulatory subunit (IKKγ). Upon its activation, the IKK complex phosphorylates IκB proteins (IκBα, IκBβ phospho Ser32, 36), which marks them for ubiquitination and degradation. This enables the NFκB complex to translocate to the nucleus where it regulates transcription by binding to DNA.

Immunogen: This NFκB p50 antibody was raised against a synthetic peptide corresponding to amino acid residues 363-376 of human NFκB p50.

Buffer: PBS containing 0.2% gelatin, 0.01% thimerosal and 0.1% sodium azide. Thimerosal and sodium azide are highly toxic.

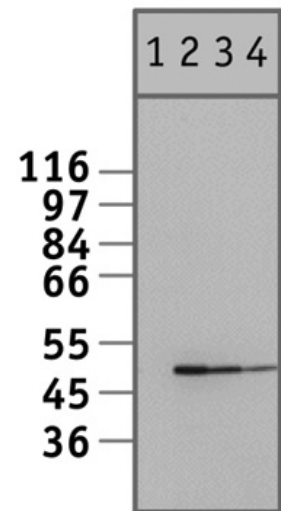
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:1,000 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.



NFκB p50 antibody tested by Western blot.

Detection of NFκB p50 by Western blot analysis. NFκB p50 can be detected in nuclear extracts derived from Raji cells using NFκB p50 rabbit polyclonal antibody. Lanes 2-4 demonstrate progressive titration of the antibody at 1:100, 1:500 and 1:1,000 dilutions, while lane 1 illustrates elimination of the signal in the presence of the immunizing peptide.