

NFκB p65 antibody (pAb)

Catalog Nos: 39283, 39284

RRID: AB_2793217

Application(s): WB

Reactivity: Human

Volumes: 200 μl, 10 μl

Purification: Affinity Purified

Host: Rabbit

Isotype: IgG

Molecular Weight: 65 kDa

Background: NFκB p65 (Nuclear factor NF-kappa-B p65 subunit also known as NFκB p65 or RelA) is a member of the Rel 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 that bind DNA and activate transcription. 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, IκBα 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 p65 antibody was raised against a peptide immunogen corresponding to N-terminal p65.

Buffer: Immunoaffinity purified antibody formulated in 30% glycerol containing 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

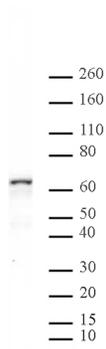
Applications Validated by Active Motif:

WB: 1:500 - 1:2,000 dilution

CUT&RUN: 1 μl per 50 μl reaction

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 p65 pAb tested by Western blot.

NFκB p65 detection by Western blot. The analysis was performed using 20 µg HeLa nuclear extract and NFκB p65 pAb at a 1:2,000 dilution.

NFκB p65 antibody (pAb) tested by CUT&RUN

CUT&RUN was performed using 500,000 HeLa nuclei and sequenced using 38 base-pair, paired-end reads on the Illumina NovaSeq. Data was collected from 16 million reads, and NFκB data is shown for Chromosome 19.

