NFkB p65 polyclonal antibody

Nuclear factor kappa-B (NFk) B) functions as a sequence specific transcriptional activator that binds to the intronic enhancer of kappa light chain gene in B lymphocytes. NF-kB is a heterodimer consisting of a 50 kDa DNA binding subunit (p50) and a 65 kDa transactivation subunit (p65/RelA) (ref. 3), both of which exhibit sequence homology to the protooncogene c-Rel. p50 has an isoform called p49/p52, and both proteins derive from the amino-terminal of precursor protein p105 and p100, respectively. The p50/p65 heterodimer remains in inactive form in the cytosol as a complex with its inhibitor, IkB. Upon stimulation of cells by a variety of stimuli, including lipopolysaccharide (LPS), pro-inflammatory cytokines like interleukin-1 and tumor necrosis factor, and viral infection, IkB becomes phosphorylated and degraded by the proteosome. The active NF-kB heterodimer translocates into the nucleus and induces gene expression.

This antibody is covered by our Worry-Free Guarantee.

Citations: 5

View Online »

Ordering Information

Order Online »

ADI-KAS-TF110-D	50µg
ADI-KAS-TF110-F	200μg

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name RELA

Application WB

Application NotesDetects a band of ~65kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

GenBank ID M62399

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of human p65

NF-κB. The sequence is completely conserved in mouse

NF-κB p65.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/ConditionsWestern Blot (1:1,000, colorimetric)Suggested

dilutions/conditions may not be available for all

applications. Optimal conditions must be determined

individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Dog, Guinea pig, Human, Monkey, Mouse,

Porcine, Rabbit, Rat

UniProt ID Q04206

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com