
NF-kB p65 Polyclonal Antibody

(Catalog # A71092)

Background

NF-kappa-B is a ubiquitous transcription factor involved in several biological processes. It is held in the cytoplasm in an inactive state by specific inhibitors. Upon degradation of the inhibitor, NF-kappa-B moves to the nucleus and activates transcription of specific genes. NF-kappa-B is composed of NFKB1 or NFKB2 bound to either REL, RELA, or RELB. The most abundant form of NF-kappa-B is NFKB1 complexed with the product of this gene, RELA. Four transcript variants encoding different isoforms have been found for this gene.

Description

NF-kB p65 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q04206

Purification

Affinity Purification

Immunogen

A synthetic peptide of human NF-kB p65

Storage

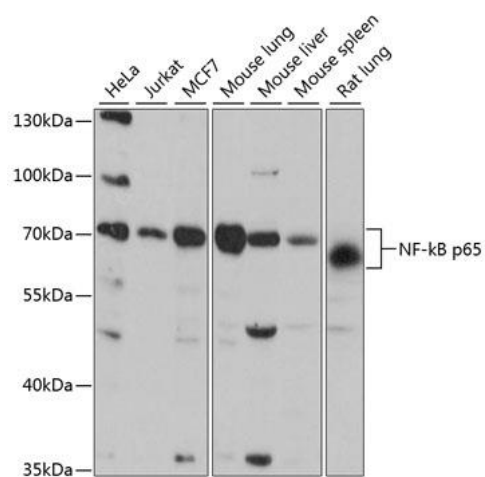
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

RELA; NFKB3; p65; transcription factor p65

Application

WB, IF; Recommended dilution: WB: 1:500-1:2000, IF:1:50-1:200



Western blot analysis of extracts of various cell lines, using NF-κB p65 Polyclonal Antibody at 1:3000 dilution.
Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 3min.