

---

## NF-kB p65 Polyclonal Antibody

(Catalog # A71094)

---

### 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 Purified

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 50-180 of human NF-kB p65/RelA (NP\_068810.3)

### Storage

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

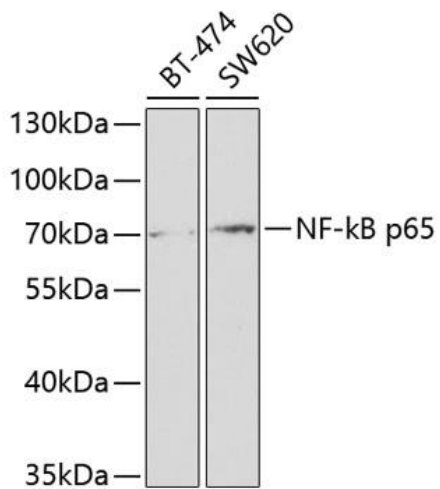
### Alternative Names

RELA; NFKB3; p65; transcription factor p65

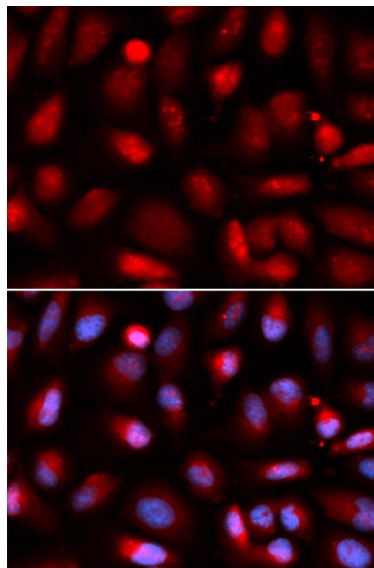
### Application

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

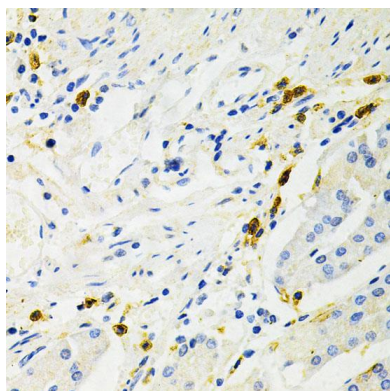
---



Western blot analysis of extracts of various cell lines, using NF-kB p65 Polyclonal Antibody at 1:1000 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.



Immunofluorescence analysis of U2OS cells using NF-kB p65 Polyclonal Antibody.



Immunohistochemistry of paraffin-embedded human stomach using NF-kB p65 Polyclonal Antibody at dilution of 1:200 (40x lens).