

## Anti-Rel A

Rabbit Polyclonal Antibody

**Catalog # R29-63BR**

Lot # M132-6

### Cited Applications

- Western blot (1:1000)

*Ideal working dilutions for each application should be empirically determined by the investigator.*

### Specificity

Recognizes the Rel A (NFκB/p65) protein

### Cross Reactivity

- Western blot human and mouse Rel A

*Rel A from other species may also be detectable*

### Host

Rabbit

### Immunogen

Synthetic peptide corresponding to amino acids 527-531 of human Rel A

### Formulation

PBS, pH7.4, 0.02%NaN<sub>3</sub>, 50% glycerol

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

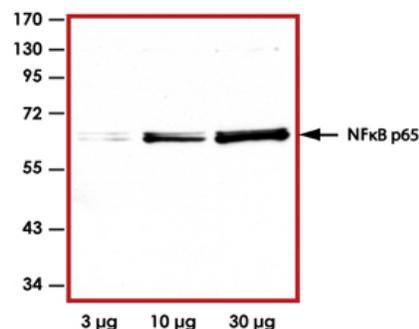
### Scientific Background

Transcription factors of the nuclear factor κB (NF-κB)/Rel family play a pivotal role in inflammatory and immune responses in unstimulated cells, NF-κB is sequestered in the cytoplasm by its inhibitory proteins, the IκBs (1). NF-κB-activating agents can induce the phosphorylation of IκB's, targeting them for rapid degradation through an ubiquitin-proteasome pathway, releasing NF-κB to enter the nucleus, where it regulates gene expression.

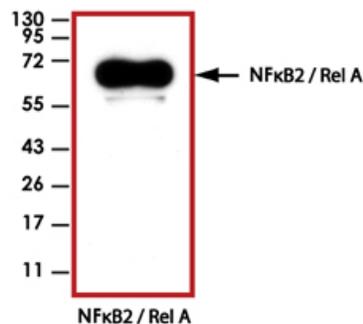
### References

1. Baeuerle, P. et al: Function and activation of NF-kappa B in the immune system. *Annu Rev Immunol.* 1994; 12: 141-79.

### Sample Data



Representative Western blot with Anti-Rel A (1:1000) using 3 µg, 10 µg, and 30 µg of 3T6 cell lysate.



Representative Western blot with Anti-Rel A (1:1000) using 20 ng human recombinant Rel A.

## Anti-Rel A

Rabbit Polyclonal Antibody

Catalog Number

R29-63BR

Specific Lot Number

M132-6

Purification

Affinity chromatography

Concentration

1.0 µg/µL

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**