

# NFκB p65 polyclonal antibody

Nuclear factor kappa-B (NFκ) B) functions as a sequence specific transcriptional activator that binds to the intronic enhancer of kappa light chain gene in B lymphocytes. NF-κB 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, IκB. 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, IκB becomes phosphorylated and degraded by the proteasome. The active NF-κB heterodimer translocates into the nucleus and induces gene expression.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 6

[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 Notes	Detects 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/Conditions	Western 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 <a href="#">Worry-Free Guarantee</a> .



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)