

Anti-Phospho-IKKε

Rabbit Polyclonal Antibody

Catalog # I04-65R

Lot # B3216-53

Cited Applications

ELISA, WB

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

Specificity

Recognizes the human IKKε protein phosphorylated at Threonine 501

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

The antibody was produced against synthesized peptide corresponding to a region of the protein surrounding phosphorylated Thr501 conjugated to KLH

Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) Sodium Azide

Stability

1yr at -20°C from date of shipment

Scientific Background

Nuclear Factor kappa B (NF-κB) is a ubiquitous transcription factor and an essential mediator of gene expression during the activation of immune and inflammatory responses. NF-κB mediates the expression of a great variety of genes in response to extracellular stimuli. NF-κB is associated with IκB proteins in the cytoplasm of the cell, which inhibit NF-κB activity. IκB proteins are phosphorylated by an IκB kinase complex consisting of at least three proteins, IKKα, IKKβ, and IKKγ. Isolated from a cDNA library of LPS-stimulated mouse macrophage cells, a novel molecule in the IKK complex has been recently identified and designated IKKi and/or IKKe. IKKe is required for the activation of NF-κB by mitogens and T cell receptors but not by TNFα or IL-1. LPS increases the IKKe mRNA level in mouse macrophage cell lines. This protein has significant sequence homology with kinase domains of IKKα and IKKβ. Overexpression of wild type IKKe in cells phosphorylates Ser32 and Ser36 of IκBα. Anti-IKKe pT501 antibody is ideal for investigators involved in NFκappaB and apoptosis research.

References:

1. Sweeney SE, et al.: Antiviral gene expression in rheumatoid arthritis: role of IKKepsilon and interferon regulatory factor 3. *Arthritis Rheum.* 2007; 56(3): 743-752.

Anti-Phospho-IKKε

Rabbit Polyclonal Antibody

Catalog # I04-65R

Lot # B3216-53

Purification Immunoaffinity chromatography and cross adsorption

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
www.signalchem.com

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