

Anti-IKK alpha

Rabbit Polyclonal Antibody

Catalog # C51-63BR

Lot # O2109-9

Cited Applications

WB, ELISA, ICC, IF

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

Specificity

Recognizes the IKK alpha protein and has no cross response to IKK beta or IKK gamma

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

IKK alpha antibody was raised against a peptide corresponding to amino acids 699 to 715 of human IKK alpha, which differs from corresponding murine sequence by one amino acid.

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

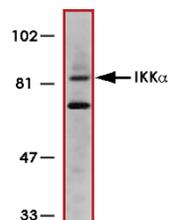
Scientific Background

IKK alpha is a serine/threonine protein kinase that phosphorylates the I-kappa-B protein which is an inhibitor of the transcription factor NF-kappa-B complex. Phosphorylation of I-kappa-B protein triggers the degradation of the inhibitor via the ubiquitination pathway, thereby activating NF-kappa-B complex (1). IKK alpha is an essential regulator of NF-kappa-B-dependent gene expression through control of promoter-associated histone phosphorylation after cytokine exposure (2). IKK alpha is a critical component of the cytoplasmic transductional-transcriptional processor leading to induction of IFN alpha production. IKK alpha is also involved in the epidermis where it antagonizes mitogenic and angiogenic signals and represses tumor progression and metastases.

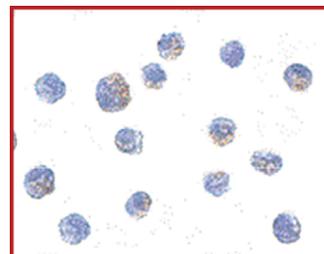
References

1. Anest, V.et.al: A nucleosomal function for I-kappa-B kinase-alpha in NF-kappa-B-dependent gene expression. Nature 423: 659-663, 2003.
2. Hoshino, K. et.al: I-kappa-B kinase-alpha is critical for interferon-alpha production induced by Toll-like receptors 7 and 9. Nature 440: 949-953, 2006.

Sample Data



Western blot analysis of IKK alpha in HeLa whole cell lysate with IKK alpha antibody at 1:500 dilution.



Immunocytochemistry of IKK alpha in HeLa cells with IKK alpha antibody at 10 µg/ml.

Anti-IKK alpha

Rabbit Polyclonal Antibody

Catalog Number

C51-63BR

Specific Lot Number

O2109-9

Purification

Affinity chromatography

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.