

Anti-IKK gamma

Rabbit Polyclonal Antibody

Catalog # I25-63R

Lot # O2109-16

Cited Applications

WB, ELISA, ICC

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

Specificity

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

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

IKK gamma antibody was raised against a peptide corresponding to amino acids 400 to 416 of human IKK gamma, which are identical to those of mouse homologue.

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

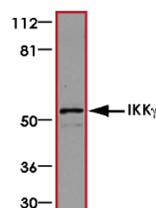
Scientific Background

IKK gamma 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. IKK gamma forms dimers and trimers and interacts preferentially with IKK beta but not IKK alpha (1). IKK gamma associates with activated ATM after the induction of DNA double-strand breaks (2). ATM phosphorylates serine-85 of IKK gamma to promote its ubiquitin-dependent nuclear export.

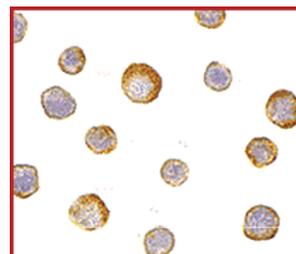
References

1. Rothwarf, D. M. et al: IKK-gamma is an essential regulatory subunit of the I-kappa-B kinase complex. *Nature* 395: 297-300, 1998.
2. Wu, Z-H. et al: Molecular linkage between the kinase ATM and NF-kappaB signaling in response to genotoxic stimuli. *Science* 311: 1141-1146, 2006.

Sample Data



Western blot analysis of IKK gamma in HeLa whole cell lysate with IKK gamma antibody at 1 µg/ml.



Immunocytochemistry of IKK gamma in HeLa cells with IKK gamma antibody at 5 µg/ml.

Anti-IKK gamma

Rabbit Polyclonal Antibody

Catalog Number

I25-63R

Specific Lot Number

O2109-16

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.