

Anti-TBK1

Rabbit Polyclonal Antibody

Catalog # T02-63R

Lot # O2109-17

Cited Applications

WB, ELISA, ICC

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

Specificity

Recognizes the TBK1 protein

Cross Reactivity

Human and Mouse

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

TBK1 antibody was raised against a synthetic peptide corresponding to 17 amino acids from near the carboxy terminus of human TBK1.

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

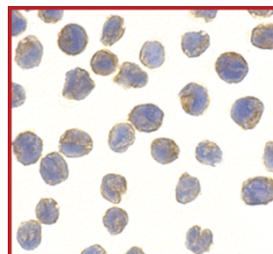
Scientific Background

TBK1, also known as NAK or NFκB-activating kinase, is an upstream protein kinase that can phosphorylate and activate the IκB kinases (1). Activation of IκB kinases allows the phosphorylation of IκB protein which is then degraded via the ubiquitination pathway. This mechanism allows the activation of the NFκB transcriptional complex. TBK1 is a specific upstream regulator of IκB kinases and can also interact and the IκB protein TANK. TBK1 is a component of the virus-activated kinase that phosphorylate IRF3 and IRF7 allowing their dimerization and translocation to the nucleus, where they induce transcription of interferon (2).

References

1. Tojima, Y. et al: NAK is an I-kappa-B kinase-activating kinase. Nature 404: 778-782, 2000.
2. Sharma, S. et al: Triggering the interferon antiviral response through an IKK-related pathway. Science 300: 1148-1151, 2003.

Sample Data



Immunocytochemistry of TBK1 in MOLT4 cells with NAK antibody at 10 µg/ml.

Anti-TBK1

Rabbit Polyclonal Antibody

Catalog Number

T02-63R

Specific Lot Number

O2109-17

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.