

IKKγ Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # I25-30G

Lot # N293-1

Product Description

Recombinant full-length human IKKγ was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_003639](#).

Gene Aliases

IKBKγ; AMCBX1; FIP-3; FIP3; Fip3p; IKK-gamma; IP; IP1; IP2; IPD2; NEMO

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

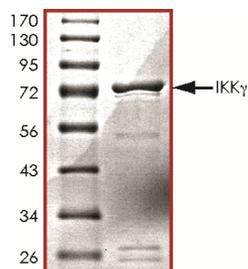
Scientific Background

IKKγ 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. IKKγ forms dimers and trimers and interacts preferentially with IKKβ but not IKKα (1). IKKγ associates with activated ATM after the induction of DNA double-strand breaks (2). ATM phosphorylates serine-85 of IKKγ 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.

Purity



The purity of IKKγ was determined to be **>95%** by densitometry. Approx. MW **~73kDa**.

IKKγ Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	I25-30G
Specific Lot Number	N293-1
Purity	>95%
Concentration	0.2 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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.