

Recombinant IKK β protein

Catalog No: 31176

Expressed In: Baculovirus

Quantity: 10 μ g

Concentration: 0.695 μ g/ μ l

Source: Human

Buffer Contents: 10 μ g of Recombinant IKK β protein in 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 20% glycerol. Protein is supplied at 0.695 μ g/ μ l.

Background: IKK β (I κ B Kinase beta subunit) – The I κ B Kinase (IKK) complex is the key enzyme involved in the activation and translocation of NF κ B (NF κ B p50 and NF κ B p65). The IKK complex is composed of two catalytic subunits (IKK α and IKK β) and a regulatory subunit (IKK γ).

NF κ B signaling is negatively regulated by the sequestration of the NF κ B complex in the cytoplasm by its association with the I κ B family of inhibitory proteins. Upon its activation, the IKK complex phosphorylates I κ B α at Ser32 and Ser36. This results ultimately in the degradation of I κ B and the activation and nuclear translocation of NF κ B.

While both IKK-alpha and IKK-beta subunits are Ser/Thr kinases, IKK beta appears to be the principal kinase, as IKK α is not required for activation of the IKK complex and the subsequent degradation of I κ B α . Once activated, IKK β autophosphorylates, which decreases IKK activity, thereby preventing a prolonged activation of inflammation.

Protein Details: Recombinant human IKK β protein was produced using baculovirus infected Sf9 cells. The protein was made against amino acids M1-S756, accession number NM_001556 and N-terminally fused to GST-HIS₆-Thrombin cleavage site. Purified by Immobilized Metal Affinity Chromatography.

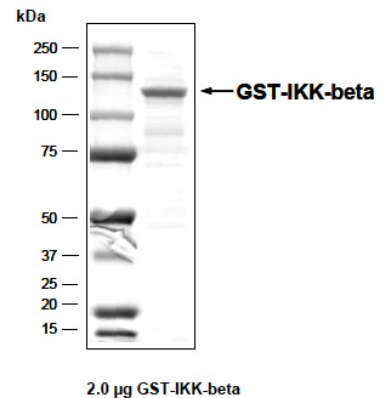
Application Notes: Recombinant IKK β is suitable for kinase assays and Western blot. The molecular weight of the protein is 120.201 kDa. The activity of the protein is 2 pmol/ μ g min.

Recommended kinase reaction conditions: 60 mM HEPES-NaOH, pH 7.5, 3 mM MgCl₂, 3 mM MnCl₂, 3 μ M Na-orthovanadate, 1.2 mM DTT, 50 μ g/ml PEG_{20,000}, ATP (variable), Substrate: I κ B α derived peptide (R11-DDRHDSGLDSMKD) at 40 μ g/ml, Recombinant IKK β at 2 μ g/ml.

Kinase activity may vary depending on the substrate and reaction conditions used.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

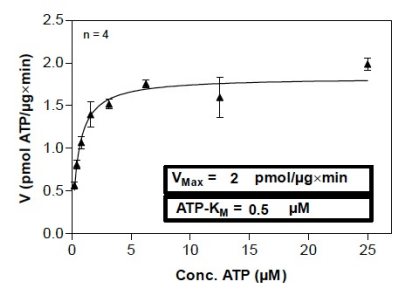
This product is for research use only and is not for use in diagnostic procedures.



Recombinant IKK β protein Coomassie Stain

Coomassie stained gel of 2 μ g of IKK β protein.

Determination of K_m value for ATP:



IKK β activity assay.

Recombinant IKK β activity measured using a kinase assay.