

Recombinant IRAK-4 protein

Catalog No: 31179

Expressed In: Baculovirus

Quantity: 10 µg

Source: Human

Buffer Contents: 10 µg of Recombinant IRAK-4 protein in 50 mM Tris-HCl, pH 8.0; 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol. Protein is supplied at 0.195 µg/µl.

Protein Details: Recombinant human IRAK-4 protein was produced using baculovirus infected Sf9 cells. The protein was made against amino acids A104-S460, accession number AF445802 and N-terminally fused to GST-HIS6-Thrombin cleavage site. Purified by GSH-agarose affinity purification.

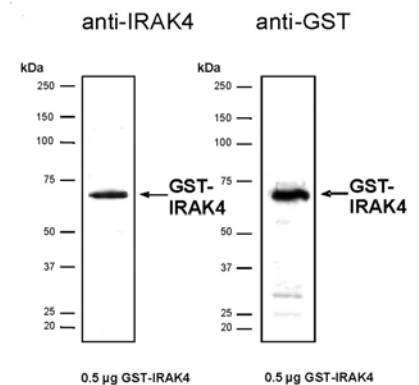
Application Notes: Recombinant IRAK-4 is suitable for kinase assays and Western blot. The molecular weight of the protein is ~69.898 kDa. The activity of the protein is ~ 574 pmol/µg min.

Recommended kinase reaction conditions: 60 mM HEPES-NaOH, pH 7.5, 3 mM MgCl₂, 3 mM MnCl₂, 3 µM Na-orthovanadate, 1.2 mM DTT, ATP (variable), 2.5 µg/50 µl PEG20.000, Substrate: Histone H2B 5 µg/50 µl, Recombinant IRAK-4 : 20 ng/50 µl.

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

Storage and Guarantee: 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 IRAK-4 protein tested by SDS-PAGE gel.

0.5 µg of IRAK-4 protein was analyzed using SDS/PAGE followed by Western blot with either an anti-IRAK-4 or anti-GST antibody.