

Recombinant FGF-R3 protein

Catalog No: 31170

Lot No: 02606001

Expressed In: Baculovirus

Quantity: 10 µg

Concentration: 0.103 µg/µl

Source: Human

Buffer Contents: 10 µg of Recombinant FGF-R3 protein in 50 mM Tris-HCl, pH 8.0; 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol.

Protein Details: Recombinant human FGF-R3 protein was produced using baculovirus infected Sf9 cells. The protein was made against amino acids E322-T806, accession number NM_000142 and N-terminally fused to a GST-HIS6-Thrombin cleavage site. Purified by GSH-agarose affinity purification.

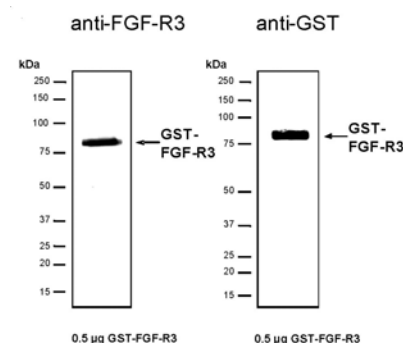
Application Notes: Recombinant FGF-R3 is suitable for kinase assays and Western blot. The molecular weight of the protein is ~82.423 kDa. The activity of the protein is ~8 pmol/µg min.

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: PolyEY 0.5 µg/50 µl, Recombinant FGF-R3: 100 ng/50 µl.

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 FGF-R3 protein tested by SDS-PAGE gel.

0.5 µg of FGF-R3 protein was analyzed using SDS/PAGE followed by Western blot with either an anti-FGF-R3 or anti-GST antibody.