

Recombinant B-RAF protein

Catalog No: 31161

Lot No: 02606001

Expressed In: Baculovirus

Quantity: 10 µg

Concentration: 0.294 µg/µl

Source: Human

Buffer Contents: 10 µg of Recombinant B-RAF protein in 50 mM Tris-HCl, pH 8.0; 100 mM NaCl, 5 mM DTT, 4 mM reduced glutathione, 20% glycerol. Protein is supplied at 0.294 µg/µl.

Protein Details: Recombinant human B-RAF protein was produced using baculovirus infected Sf9 cells. The protein was against amino acids Gln417-His766, accession number NM_004333.2 and N-terminally fused to a GST-HIS6-Thrombin cleavage site. Purified by GSH-agarose affinity purification.

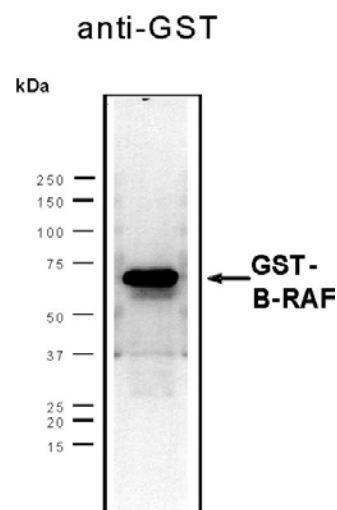
Application Notes: Recombinant B-RAF is suitable for kinase assays and Western blot. The molecular weight of the protein is ~69.127 kDa. The activity of the protein is ~ 78 pmol/µg min.

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: MEK1-DN, 5 µg/50 µl, Recombinant B-RAF: 25 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.



500 ng GST-B-RAF

Recombinant B-RAF protein tested by SDS-PAGE gel.

0.5 µg of B-RAF protein was analyzed using SDS/PAGE followed by Western blot with an anti-GST antibody.