

Recombinant JAK2 protein

Catalog No: 31181

Expressed In: Baculovirus

Quantity: 10 µg

Concentration: 0.091 µg/µl

Source: Mouse

Buffer Contents: 10 µg of Recombinant JAK2 protein in 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 4 mM reduced glutathione, 20% glycerol. Protein is supplied at 0.091 µg/µl.

Protein Details: Recombinant mouse JAK2 protein was produced using baculovirus infected Sf9 cells. The protein is a C-terminal fragment containing amino acids P717-G1132, accession number NP_004963.1 fused to an N-terminal GST-HIS₆ tag with a Thrombin cleavage site. Purified by GST-affinity chromatography.

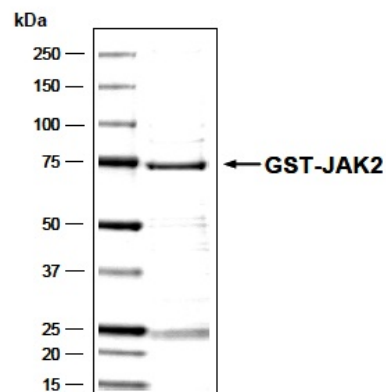
Application Notes: Recombinant JAK2 is suitable for kinase assays and Western blot. The molecular weight of the protein is 78.512 kDa. The activity of the protein is 24 pmol/µg min. ATP-K_M: 2.1 µM

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), Substrate: TRK-C-derived Peptide (R11-VYSTDYRFLNPS) at 50 µg/ml, Recombinant JAK2 at 4 µg/ml.

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.

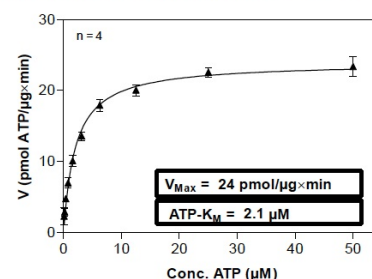
This product is for research use only and is not for use in diagnostic procedures.



Recombinant JAK2 protein Coomassie Stain

Coomassie stained gel of 2 µg of JAK2 protein.

Determination of V_{max} and K_M value for ATP



JAK2 activity assay.

Recombinant JAK2 activity measured using a kinase assay.