

## Recombinant JAK3 protein

**Catalog No:** 31182

**Expressed In:** Baculovirus

**Quantity:** 10 µg

**Source:** Human

**Buffer Contents:** 10 µg of Recombinant JAK3 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 0.213 µg/µl.

**Protein Details:** Recombinant human JAK3 protein was produced using baculovirus infected Sf9 cells. The protein was made against amino acids I781-S1124, accession number NM\_000215 and N-terminally fused to GST-HIS6-Thrombin cleavage site. Purified by GSH-agarose affinity purification.

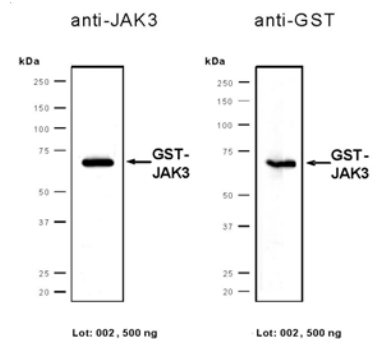
**Application Notes:** Recombinant JAK3 is suitable for kinase assays and Western blot. The molecular weight of the protein is ~68.379 kDa. The activity of the protein is ~ 8 pmol/µg min.

Recommended kinase reaction conditions: 60 mM HEPES-NaOH, pH 7.5, 3 mM MgCl<sub>2</sub>, 3 mM MnCl<sub>2</sub>, 3 µM Na-orthovanadate, 1.2 mM DTT, ATP (variable), 2.5 µg/50 µl PEG20.000, Substrate: R11-TRK-C-derived Peptide (R11-VYSTDYRLFNPS) 2.5 µg / 50 µl, Recombinant JAK3 : 200 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 JAK3 protein tested by SDS-PAGE gel.**

0.5 µg of JAK3 protein was analyzed using SDS/PAGE followed by Western blot with either an anti-JAK3 or anti-GST antibody.