

JAK2 (JH1, JH2 domain), GST-Tag, Avi-Tag, Biotin (Dephosphorylated) Recombinant

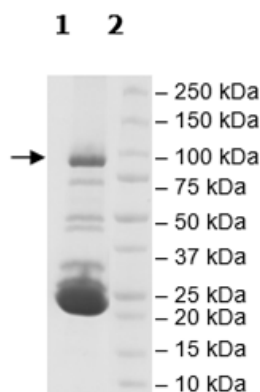
Catalog: 79495
Lot: 200623

Product Information

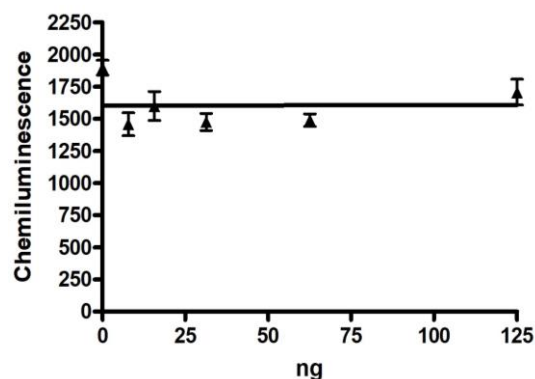
Description:	Recombinant human JAK2 containing the catalytic and the pseudo kinase domain (Just Another Kinase 2), encompassing amino acids 513-1132(end). This construct contains an N-terminal GST-Tag and a C-terminal Avi-Tag™ and was co-expressed with PTP1B. The recombinant protein was affinity purified.
Species	Human
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
Construct:	JAK2 (GST-1-1132(end)-Avi)-(Biotin)
Concentration:	0.49 mg/ml
Expression System:	HEK293
Purity:	15%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT, and 16 mM glutathione
MW:	100 kDa
Genbank Accession:	NM_004972
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	Assay was performed according to TYR Phosphorylation Assay Kit (BPS Bioscience #79502), with both phosphorylated JAK2 (standard) and JAK2 titrated from 0 ng to 500 ng.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



JAK2 Activity



Quality Control Data

Biotin-Avidin Pulldown

