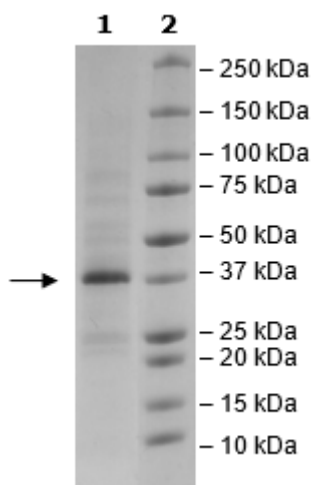


Product Information

Construct:	JAK2 (JH1 domain) (His-808-1132(end))
Concentration:	0.52 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 0.04% Tween-20 and 20% Glycerol
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_004972
MW:	43 kDa
Purity:	71%
Specific Activity:	124 pmol/min/μg
Assay Conditions:	Assay was done according to JAK2 (Janus Kinase 2) Assay Kit (BPS Bioscience #79520) with 20 μM ATP instead of 10 μM ATP and various amounts of JAK2 enzyme.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



JAK2 Activity

