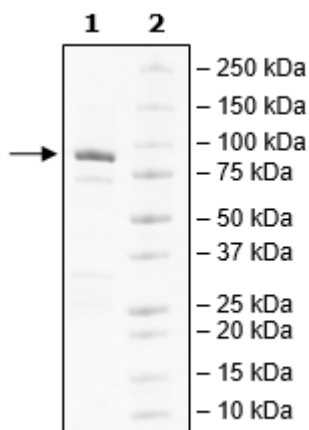


## Product Information

<b>Construct:</b>	JAK2 (JH1, JH2 Domain) (His-GST-532-1132(end))
<b>Concentration:</b>	0.30 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	45 mM Tris-HCl, pH 8.0, 300 mM NaCl, 2.4 mM KCl, 0.04% Tween-20, 3 mM DTT, 10% glycerol and 250 mM Imidazole
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_004972
<b>MW:</b>	98 kDa
<b>Purity:</b>	80%
<b>Assay Conditions:</b>	Assay was done in a kinase buffer containing 1 mM DTT using Poly-(Glu4:Tyr) substrate (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using 50 μl of Kinase-Glo reagent (Promega) per well.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



JAK2 (JH1, JH2 domain) Activity

