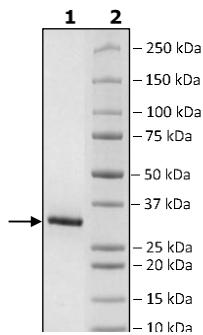


Product Information

Construct:	JAK2(W659A, W777A, F794H) (536-812-Th-His)
Mutations:	W659A, W777A, F794H
Concentration:	2.25 mg/ml
Species:	Human
Formulated In:	20 mM Tris HCL, pH 8.5, 0.5 mM TCEP and 10% 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:	33 kDa
Purity:	≥90%
Assay Conditions:	Assay was done according to the JAK2 JH2 Pseudokinase Domain Inhibitor Screening Assay Kit (BPS Bioscience #78111) with JAK2 titrated from 1000 nM to 1.5 nM.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Binding of JH2 Probe 1 to JAK2 JH2 Domain

