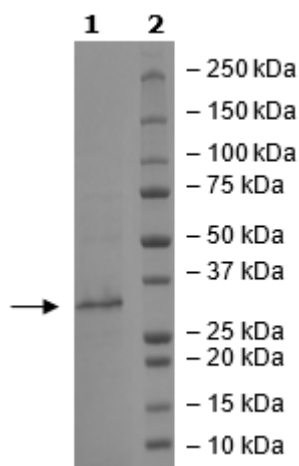


Product Information

Construct:	JAK2 (V617F, W659A, W777A, F794H) (536-812-Th-His)
Mutations:	V617F, W659A, W777A, F794H
Concentration:	1.09 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.5mM TCEP, 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:	33 kDa
Purity:	≥90%
Assay Conditions:	Assay was done according to JAK2-JH2 Pseudokinase Domain (V617F, W659A, W777A, F794H) Inhibitor Screening Assay Kit (BPS Bioscience #78209) with Jak2-JH2 titrated from 288 nM to 0 nM.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



JAK2-JH2 Activity

