

Product Information

Construct:	JAK2 (513-827-His-Avi)-(Biotin)
Concentration:	8.56 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 300 mM NaCl, 2.2 mM KCl, 20% glycerol, 0.2mM TCEP
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:	39 kDa
Purity:	>90%
Notes:	This protein is enzymatically biotinylated using Avi-tag™ technology. Biotinylation is confirmed to be ≥90%.

Quality Control Data

4-20% SDS-Page Coomassie Staining

