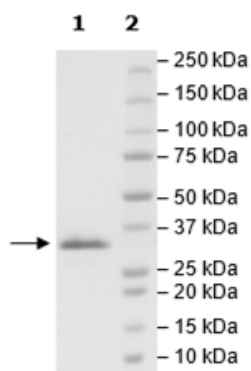


Product Information

<b>Construct:</b>	JAK2 (V617F, W659A, W777A, F794H) (536-812-His-Avi)-(Biotin)
<b>Mutations:</b>	V617F, W659A, W777A, F794H
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
<b>Concentration:</b>	0.66 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	45 mM Tris-HCl, pH 8.0, 270 mM NaCl, 2 mM KCl, 20% glycerol, 0.5 mM TCEP
<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>	35 kDa
<b>Purity:</b>	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown

