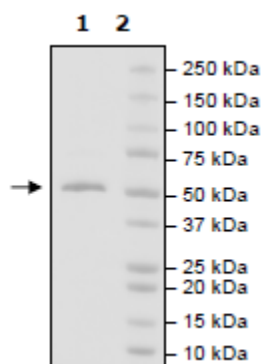


Product Information

Construct:	His-Avi(biotin)-FGFR3(397-806(end)) (dephosphorylated)
Concentration:	1.63 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT
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_000142
MW:	47 kDa
Purity:	≥90%
Notes:	This protein was co-expressed with PTP1B for dephosphorylation. Phosphorylation level confirmed to be <5%. Biotinylation is confirmed to be ≥90%.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown

