

Product Information

Construct:	FGFR3 (N540K) (His-397-806(end)) / PTP1B (GST-Th-1-435(end))-(Biotin)
Concentration:	0.17 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:	70%
Notes:	This protein was co-expressed with PTP1B for dephosphorylation. Phosphorylation level confirmed to be <5%. This protein was purified by mono-Avidin with biotin removed by buffer exchange. Biotinylation is confirmed to be ≥90%.

Quality Control Data

4-20% SDS-Page Coomassie Staining

