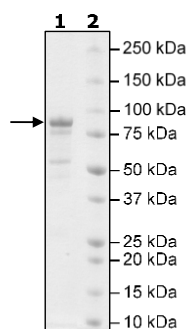


Product Information

Construct:	VEGFR3 (FLT4) (GST-798-1298)
Concentration:	0.38 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_182925
MW:	82 kDa
Purity:	63%
Assay Conditions:	Assay was done according to the VEGFR3 (FLT4) Kinase Assay Kit (BPS Bioscience #79738) with VEGFR3(FLT4) titrated from 400 ng/rxn – 0.625 ng/rxn with 20 μ M ATP instead of 10 μ M.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



VEGFR3 (FLT4) Activity

