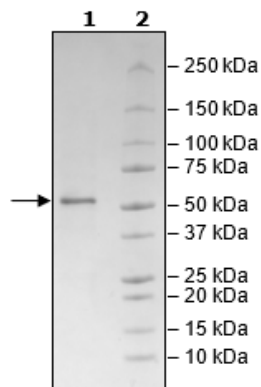


Product Information

Construct:	Flt3 (His-564-993(end))
Concentration:	0.41 mg/ml
Species:	Human
Formulated In:	25 mM Tris-HCl pH 8.0, 69 mM NaCl, 1.35 mM KCl, 0.025% Tween-20, 63 mM imidazole, 50% glycerol, and 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_004119
MW:	55 kDa
Purity:	≥90%
Assay Conditions:	Assay was done according to FLT3 Kinase Assay Kit (BPS Bioscience #79797) with 20 μM ATP and FLT3 titrated from 75 ng/rxn to 0 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



FLT3 Activity

