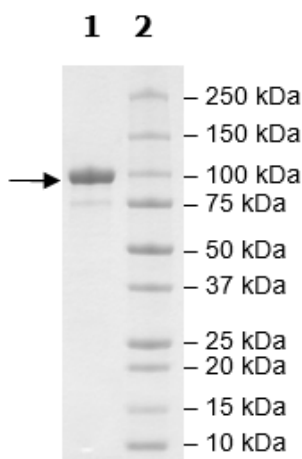


Product Information

Construct:	VEGFR2 (KDR) (GST-805-1356(end))
Concentration:	0.11 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, 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:	AF035121
MW:	88 kDa
Purity:	70%
Specific Activity:	470 pmol/min/g
Assay Conditions:	Assay was done in Kinase buffer containing 0.2 mM DTT using Poly-(Glu4:Tyr)-biotin substrate (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



VEGFR2 Activity

