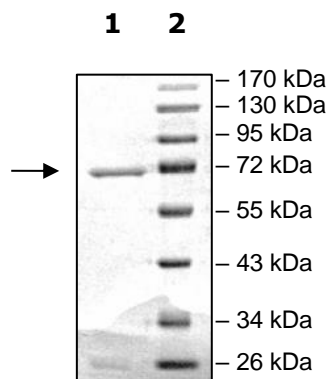


## Product Information

<b>Construct:</b>	EPHA1 (GST-569-end)
<b>Concentration:</b>	0.10 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.25 mM DTT, 0.1 mM PMSF, 25% glycerol
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_005232
<b>MW:</b>	72 kDa
<b>Purity:</b>	≥90%
<b>Specific Activity:</b>	28 pmol/min/μg
<b>Assay Conditions:</b>	Kinase activity was measured using a Poly-Glu/Tyr(4:1) substrate at 1mg/ml. The protein kinase was diluted to 0.1 μg/ μl in buffer containing 50ng/μl BSA. Increasing amounts of the protein kinase were mixed with 0.25 mg/ml substrate in 20 μl final volume. The blank was determined from a “no substrate” sample. The reaction was initiated by addition of 5 μl of [33P]-ATP diluted in kinase buffer: 6ml kinase buffer containing 1 mCi [33P]-ATP, 0.25 mM ATP, 25 mM MOPS, 12.5 mM β-glycerophosphate, 20 mM MgCl <sub>2</sub> , 12.5 mM MnCl <sub>2</sub> , 5 mM EGTA, 2 mM EDTA, 0.25 mM fresh DTT, pH 7.2. After incubating for 30°C for 15 minutes, the reaction was terminated by spotting 20 μl of the mixture onto phosphocellulose paper strips that were fixed in 1% phosphoric acid and washed three times. Radioactivity was determined using a scintillation counter.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

