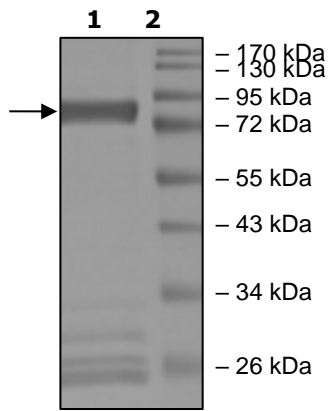


## Product Information

<b>Description:</b>	Recombinant human FGFR1 (Fibroblast Growth Factor Receptor 1), encompassing amino acids 399-822 (end) and containing the kinase domain. This construct contains an N-terminal GST-tag. This recombinant protein was affinity purified and is kinase active.
<b>Species:</b>	Human
<b>Construct:</b>	FGFR1 (GST-399-822)
<b>Concentration:</b>	0.10 mg/ml
<b>Expression System:</b>	Sf9
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<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, and 25% glycerol
<b>MW:</b>	73 kDa
<b>Genbank Accession:</b>	NM_023110
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Specific Activity:</b>	62 pmol/min/μg
<b>Assay Conditions:</b>	FGFR1 (FLT2) activity was measured by using a Poly (4:1 Glut, Tyr) Peptide Substrate, diluted in 25 mM Tris-HCl pH 7.5 to a final concentration of 1 mg/ml, in the ADP Glo™ Kinase Assay Kit (Promega #V9101). FGFR1 (FLT2) was diluted at different concentrations in 1x Kinase Assay Buffer (40 mM Tris-HCl, pH 7.4, 20 mM MgCl <sub>2</sub> , 2.5 mM MnCl <sub>2</sub> , 0.1 mg/ml BSA and 50 μM DTT). A Substrate/ATP mix was prepared by adding 1 μl of 10 mM ATP, 78 μl of 5x Kinase Assay Buffer, 80 μl of 1 mg/ml substrate and 1 μl of 1 M of MnCl <sub>2</sub> . The reaction was initiated by the addition of 3 μl of diluted FGFR1 (FLT2) to 2 μl of Substrate/ATP Mix. After a 40-minute incubation at Room Temperature (RT), the reaction was terminated by addition of ADP-Glo™ Reagent, followed by a subsequent 40-minute incubation at RT. Kinase Detection Reagent was added, and the reaction was incubated for another 30 minutes at RT. Detection of luminescence was measured using the Luminescence Module Protocol on GloMax®-Multi Microplate Multimode Reader. The Specific Activity was calculated as follows: (Corrected activity, RLU) / [(Specific activity from ADP in RLU/pmol) * (Reaction time in min) * (Enzyme amount in μg or mg)]. Corrected RLU was calculated by subtracting the blank value from all the values. The blank was determined from a “no enzyme” sample by replacing the enzyme solution with an equal volume of 1x Kinase Assay Buffer.
<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

