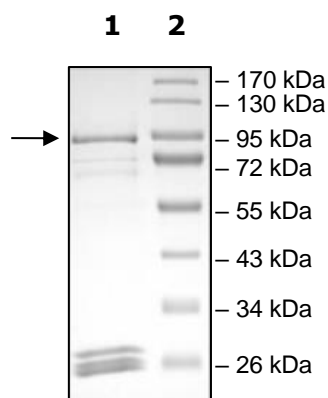


Product Information

Construct:	EGFR (GST-668-end)
Mutations:	Del746-750, T790M, C797S
Concentration:	0.1 mg/ml
Species:	Human
Formulated In:	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
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_005228
MW:	79 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	80%
Specific Activity:	17 pmol/min/μg
Assay Conditions:	Kinase activity was measured using ADP-Glo™ Kinase Assay kit (Promega; Cat# V9101) which quantifies the amount of ADP produced by phosphorylation. The ADP-Glo™ Reagent is added to terminate the reaction and quench the remaining ATP. The Kinase Detection Reagent is then added to convert ADP to ATP and to measure the newly converted ATP using a luciferase reaction. Assay: Kinase activity was measured using substrate poly-Glu,Tyr 4:1 (stock 1 mg/ml). Increasing amounts of EGFR were incubated with a final concentration of 200 μg/ml substrate in 60 mM Tris-HCl, pH 7.4, 50 mM MgCl ₂ , 4 mM MnCl ₂ , 250 μg/ml BSA, 50 μM fresh DTT, and 25 μM ATP. The total incubation time was 40 min at room temperature. The reaction was terminated by addition of an equal volume of the ADP-Glo™ Reagent supplemented with 10 mM MgCl ₂ , and the Kinase Detection Reagent was added. Phosphorylation was measured by detection of luminescence. The blank was determined from a “no kinase” sample.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

