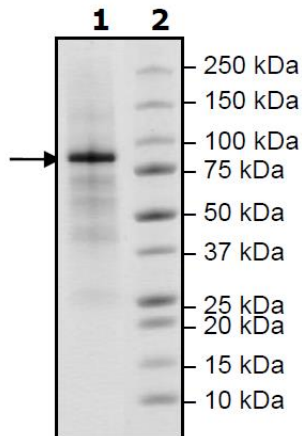


## Product Information

<b>Construct:</b>	EGFR (T790M, L858R) (GST-668-1210-His)
<b>Mutation:</b>	T790M, L858R
<b>Concentration:</b>	0.04 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 300 mM imidazole.
<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_005228
<b>MW:</b>	88 kDa
<b>Purity:</b>	70%
<b>Specific Activity:</b>	680 pmol/min/μg
<b>Assay Conditions:</b>	Assay was done in a kinase buffer containing 2 mM DTT using Poly-(Glu4:Tyr) substrate (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 40 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).
<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

