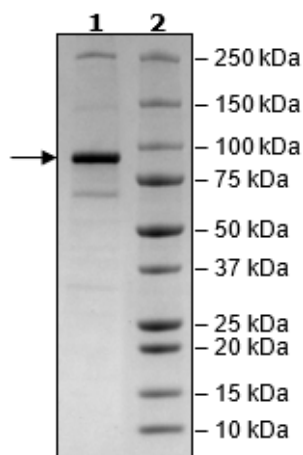


Product Information

Description:	Human recombinant EGFR (epidermal growth factor receptor, also known as HER1) encompassing amino acids 668-1210 which contains the tyrosine kinase domain. The construct contains an N-terminal GST-Tag, a C-terminal His-Tag (6xHis), and the mutation leucine to arginine at position 858. This protein was affinity purified.
Species:	Human
Construct:	EGFR (L858R) (GST-668-1210(end)-His)
Mutation:	L858R
Concentration:	0.26 mg/ml
Expression System:	Sf9
Purity:	80%
Format:	Aqueous buffer solution.
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
MW:	88 kDa
Genbank Accession:	NM_005228
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	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.
Specific Activity:	340 pmol/min/μg
Assay Conditions:	Assay was performed according to EGFR(L858R) Kinase Assay Kit (BPS Bioscience #40324) with EGFR (L858R) titrated at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



EGFR Activity

