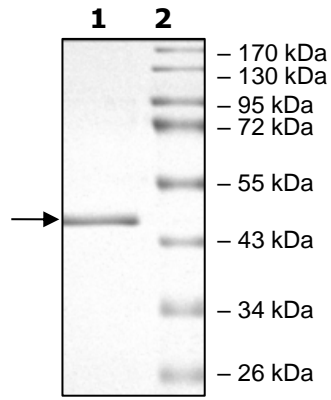


## Product Information

<b>Description:</b>	Recombinant human EGFR (epidermal growth factor receptor), encompassing amino acids (699-1022) length. This construct contains an N-terminal His-tag and mutation of interest (V948R). The recombinant protein was affinity purified and is kinase active.
<b>Species:</b>	Human
<b>Construct:</b>	EGFR (V948R) (His-669-1022)
<b>Mutation:</b>	V948R
<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 Sodium Phosphate, pH 7.0, 300 mM NaCl, 150 mM Imidazole, 0.25 mM DTT, 25% glycerol
<b>MW:</b>	48 kDa
<b>Genbank Accession:</b>	NM_005228
<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>	38 pmol/min/μg
<b>Assay Conditions:</b>	EGFR (V948R) activity was measured by using Poly (4:1 Glu, Tyr) synthetic peptide substrate diluted in 25 mM Tris-HCl buffer (pH 7.5) to a final concentration of 1 mg/ml in the ADP Glo™ assay (Promega #V9101). Reaction was initiated by mixing increasing amounts of the EGFR (V948R) with 25 μM ATP in 40 mM Tris-HCl, pH 7.4, 20 mM MgCl <sub>2</sub> , 0.1 mg/ml BSA prepared with 250 μM DTT and the 20 μg/ml substrate. After a 40-minute incubation at 37°C, the reaction was terminated by addition of the AMP-Glo™ Reagent followed by a subsequent 40-minute incubation at room temperature. Kinase Detection Reagent was then added and incubated for another 30 minutes. Detection of luminescence was measured using the Luminescence Module Protocol on GloMax®-Multi Microplate reader. The corrected activity (RLU) was calculated by removing the blank value for each sample divided by the (specific activity of ADP in RLU/pmol)*(Reaction time in min)*(Enzyme amount in μg or mg). The blank was determined from a “no kinase” sample by replacing the enzyme working solution with an equal volume of Kinase Dilution Buffer X (1X).
<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

