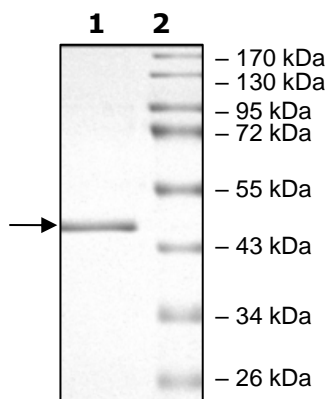


Product Information

| | |
|------------------------------|--|
| Description: | 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. |
| Species: | Human |
| Construct: | EGFR (V948R) (His-669-1022) |
| Mutation: | V948R |
| Concentration: | 0.10 mg/ml |
| Expression System: | Sf9 |
| Purity: | 90% |
| Format: | Aqueous buffer solution. |
| Formulated In: | 50 mM Sodium Phosphate, pH 7.0, 300 mM NaCl, 150 mM Imidazole, 0.25 mM DTT, 25% glycerol |
| MW: | 48 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: | 38 pmol/min/μg |
| Assay Conditions: | <p>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 MgCl₂, 0.1 mg/ml BSA prepared with 250 μM DTT and substrate with a final concentration of 50 μg/ml.</p> <p>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).</p> |
| Applications: | Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling. |

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Specific Activity

