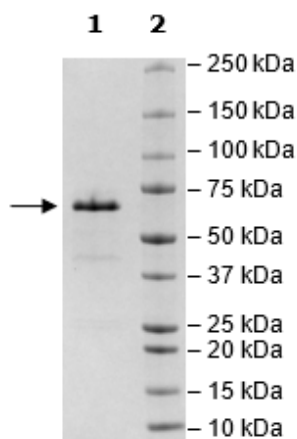


Product Information

Construct:	EGFR (Mouse) (670-1210(end)-FLAG)
Concentration:	0.13 mg/ml
Species:	Mouse
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
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_207655
MW:	62 kDa
Purity:	≥90%
Specific Activity:	≥49 pmole/min/μg
Assay Conditions:	Assay was performed according to EGFR Kinase Assay Kit (BPS Bioscience #40321).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



EGFR (Mouse) Activity

