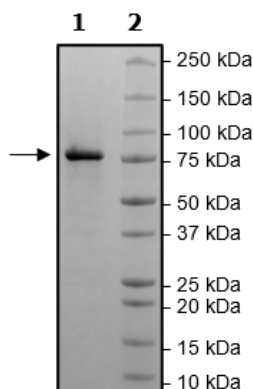


Product Information

Description:	Recombinant human CSF1R (colony stimulating factor 1 receptor), encompassing amino acids 538-972 (end) which contains the tyrosine kinase domain. This construct contains an N-terminal GST-Tag followed by a His-tag (6xHis). The recombinant protein was affinity purified.
Species	Human
Construct:	CSF1R (GST-His-Th-538-972(end))
Concentration:	1.47 mg/ml
Expression System:	Sf9
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol
MW:	79 kDa
Genbank Accession:	NM_005211
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:	≥240 pmol/min/μg
Assay Conditions:	Assay was done using ADP-Glo (Promega) reagents with 20 μM ATP, 0.2 mg/mL Poly-(Glu,Tyr) in 25 μL using Kinase Assay Buffer 1 (BPS Bioscience #79334) Reaction was performed at 30°C for 45 min.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CSF1R Activity

