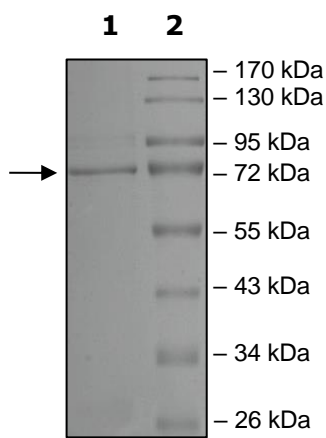


Product Information

Construct:	BTK(C481S)(His-2-end)
Mutation:	C481S
Concentration:	0.10 mg/ml
Species:	Human
Formulated In:	50 mM sodium phosphate, pH 7.0, 300 mM NaCl, 150 mM imidazole, 0.1 mM PMSF, 0.25 DTT mM, 25% glycerol.
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_000061
MW:	75 kDa
Purity:	80%
Specific Activity:	62 pmol/min/μg
Assay Conditions:	The kinase assay is performed using the ADP-Glo™ Kinase Assay kit (Promega; Cat# V9101) which quantifies the amount of ADP produced. The ADP-Glo™ Reagent is added to terminate the reaction and to deplete the remaining ATP. The Kinase Detection Reagent is then added to convert ADP to ATP and to measure the newly synthesized ATP using a luciferase reaction. Assay: kinase activity was measured using a Poly-Glu/Tyr(4:1) substrate. The reaction was initiated by mixing increasing amounts of the protein kinase with 25 μM ATP in 40 mM Tris-HCl, pH 7.4, 20 mM MgCl ₂ , 0.1 mg/ml BSA, 250 μM DTT, and 0.2 mg/ml Poly-Glu/Tyr. The reaction was terminated by addition of the ADP-Glo™ Reagent, and the Kinase Detection Reagent was added. Phosphorylation was measured by detection of luminescence. The blank was determined from a “no kinase” sample Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.
Applications:	

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

