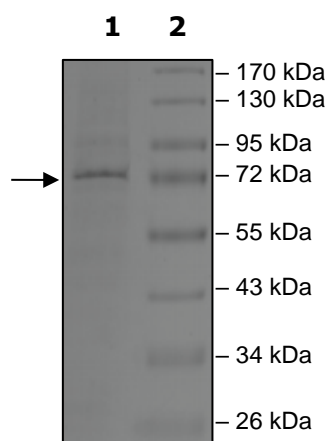


Product Information

Construct:	BTK(P190K)(His-2-658(end))
Mutation:	P190K
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, 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:	85%
Specific Activity:	220 pmol/min/μg
Assay Conditions:	Kinase activity was measured using a Poly-Glu/Tyr(4:1) substrate. Increasing amounts of the protein kinase were mixed with 0.25 mg/ml substrate in 20 μl final volume. The blank was determined from a "no substrate" sample. The reaction was initiated by addition of 5 μl of [33P]-ATP buffer containing 10 mM ATP, 1mCi [33P]-ATP, 25 mM MOPS, 12.5 mM β-glycero-phosphate, 25 mM MgCl ₂ , 2 mN EDTA, 0.25 mM fresh DTT, pH 7.2. After incubating for 30°C for 15 minutes, the reaction was terminated by spotting 20 μl of the mixture onto phosphocellulose paper strips that were fixed in 1% phosphoric acid. Radioactivity was determined using a scintillation counter.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

