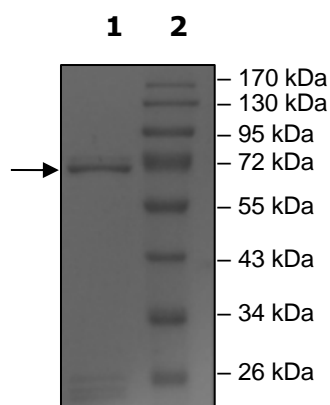


## Product Information

<b>Construct:</b>	LTK (GST-498-796)
<b>Concentration:</b>	0.10 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.25 mM DTT, 0.1 PMSF, 25% glycerol
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_002344
<b>MW:</b>	70 kDa
<b>Purity:</b>	≥90%
<b>Specific Activity:</b>	15 pmol/min/μg
<b>Assay Conditions:</b>	The activity of the complex was measured using peptide sequence (EAIYAAPFAKKK), based on the C-terminus of Abl, as a substrate (stock solution 1 mg/ml dissolved in distilled water). Assay: Increasing amounts of kinase were mixed with a final concentration of 200 μg/ml peptide substrate in a buffer consisting of 15 mM MOPS pH 7.2, 15 mM MgCl <sub>2</sub> , 7.5 mM β-glycerol-phosphate, 3 mM EGTA, 1.2 mM EDTA, 20 μg/ml BSA (bovine serum albumin) and 150 μM fresh DTT (final concentrations). The reaction was initiated by adding a [33P]-ATP (1 μCi/sample) mixed with non-radioactive ATP to reach a final concentration of 50 μM. The blank was determined from a “no substrate” sample. The reaction was incubated for 15min at 30°C and terminated by spotting 20 μl of the mixture onto phosphocellulose paper strips that were fixed in 1% phosphoric acid and washed three times. Radioactivity was determined using a scintillation counter.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

## 4-20% SDS-Page Coomassie Staining



## Specific Activity

