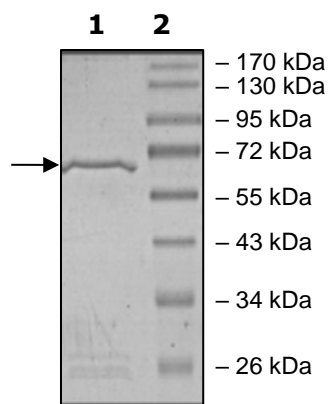


## Product Information

<b>Construct:</b>	WNK1 (GST-181-507)
<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 mM 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_018979
<b>MW:</b>	67 kDa
<b>Purity:</b>	90%
<b>Specific Activity:</b>	25 pmol/min/μg
<b>Assay Conditions:</b>	Kinase activity was measured using Myelin Basic Protein (MBP) substrate diluted in water to a final concentration of 1 mg/ml. Increasing amounts of kinase were mixed with a final concentration of 200 μg/ml peptide substrate in a buffer consisting of 5 mM MOPS pH 7.2, 5 mM MgCl <sub>2</sub> , 2.5 mM β-glycerol-phosphate, 1 mM EGTA, 0.4 mM EDTA, 50 ng/μl BSA (bovine serum albumin) and 0.25 mM fresh DTT (final concentrations). The reaction was initiated by adding 5 μl [33P]-ATP (1 μCi/sample) mixture containing 0.25 mM non-radioactive ATP. The blank was determined from a “no substrate” sample. The reaction was incubated for 15 minutes 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

