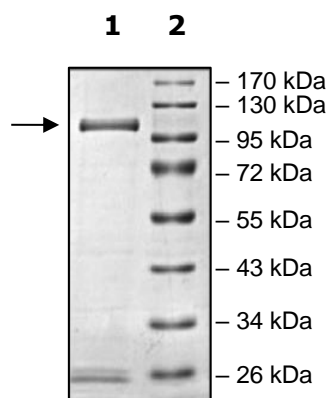


Product Information

Construct:	MSSK1 (GST-Full Length)
Concentration:	0.1 mg/ml
Species:	Human
Formulated In:	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
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	BC117124
MW:	98 kDa
Purity:	80%
Specific Activity:	210 pmol/min/μg
Assay Conditions:	Kinase activity was measured using PKCtide sequence ERMRPKRQGSVRRRV (stock solution 1 mg/ml dissolved in distilled water). 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 ₂ , 2.5 mM β-glycerol-phosphate, 1 mM EGTA, 0.4 mM EDTA, 50 μg/ml BSA (bovine serum albumin) and 50 μM fresh DTT (final concentrations). The reaction was initiated by adding [33P]-ATP (1 μCi/sample) mixed with non-radioactive ATP (final concentration of 10 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.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

