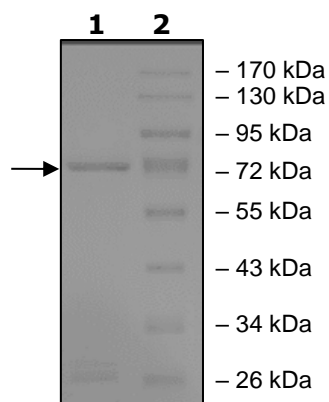


Product Information

Construct:	MEK1 (GST-Full length)
Concentration:	0.10 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:	NM_002755
MW:	72 kDa
Purity:	85%
Specific Activity:	280 pmol/min/μg
Assay Conditions:	<p>Kinase activity was measured by first activating ERK1 at 0.2 μg/μl with MEK1. Next, Myelin Basic Protein (MBP) was added as the substrate diluted in water to a final concentration of 1 mg/ml.</p> <p>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 MgCl₂, 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 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.</p> <p>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.</p>
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

