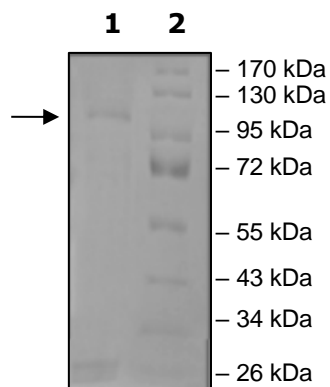


Product Information

Construct:	MEKK3 (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:	NM_002401
MW:	105 kDa
Purity:	75%
Specific Activity:	23 pmol/min/μg
Assay Conditions:	MEKK3 kinase activity was measured using MBP (Myelin Basic Protein) as substrate. Increasing amounts of the protein kinase were mixed with a final concentration of 0.25 mg/ml MBP in a buffer consisting of 5 mM MOPS pH 7.2, 2.5 mM β-glycerol-phosphate, 5 mM MgCl ₂ , 1 mM EGTA, 0.4 mM EDTA, 25 μg/ml BSA (bovine serum albumin), and 100 μM DTT (final concentrations). The reaction was initiated by adding [33P]-ATP (1 μCi/sample) mixed with non-radioactive ATP (final concentration of 50 μM). 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. The blank was determined from a “no substrate” sample.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

