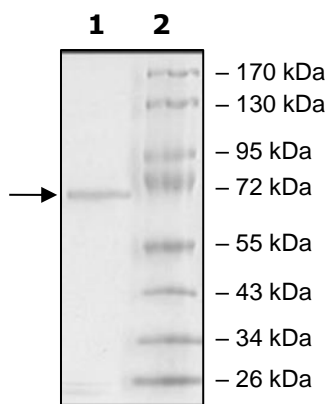


Product Information

**Construct:** p38 $\gamma$  (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\_002969  
**MW:** 71 kDa  
**Purity:** 90%  
**Specific Activity:** 220 pmol/min/ $\mu$ g  
**Assay Conditions:** p38 $\gamma$  kinase activity was measured P38 Sub synthetic peptide (IPTTPITTTYFFFKKK) diluted in water to a final concentration of 1 mg/ml. Increasing amounts of kinase were mixed with a final concentration of 200  $\mu$ g/ml peptide substrate in a buffer consisting of 5 mM MOPS pH 7.2, 2.5 mM  $\beta$ -glycerol-phosphate, 1 mM EGTA, 5 mM MgCl<sub>2</sub>, 1 mM EGTA and 0.4 mM EDTA with 50 ng/ $\mu$ l BSA (bovine serum albumin) and .25 mM fresh DTT (final concentrations). The reaction was initiated by adding 5  $\mu$ l [33P]-ATP (1  $\mu$ Ci/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  $\mu$ l 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

