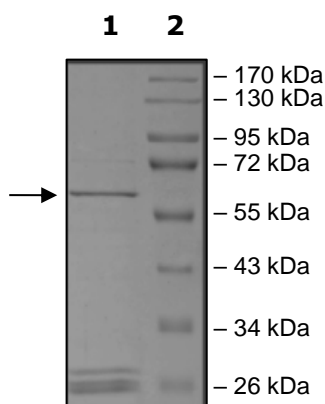


Product Information

Construct:	PP1A (GST-Full Length)
Concentration:	0.05 mg/ml
Species:	Human
Formulated In:	20 mM MOPS, pH 7.5, 300 mM NaCl, 10 mM glutathione, 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_002708
MW:	62 kDa
Purity:	85%
Specific Activity:	264 pmol/min/μg
Assay Conditions:	<p>Kinase activity was measured using BIOMOL® Green (Enzo; Cat# AK111) a colorimetric assay that quantifies the amount of free phosphates produced by the PP1A reaction. A 1:1 serial dilution of phosphate standard was prepared in a 1:4 freshly diluted Phosphatase Dilution Buffer II buffer containing .2% 2-mercaptoethanol and 65 ng/μl BSA to a final volume of 25 μl. A blank of diluted Phosphate Dilution Buffer II without phosphate was included.</p> <p>Assay: Kinase activity was measured using substrate thr-phosphopeptide synthetic substrate (KRT(p)IRR) diluted in distilled water with a concentration of 1 mg/ml. Increasing amounts of PP1A kinase were incubated with the KRT(p)IRR substrate with a final concentration of 0.2 mg/ml in Phosphatase Dilution Buffer II buffer containing 0.2% 2-mercaptoethanol and 65 ng/μl BSA in a final volume of 25 μl. The reaction was then incubated at 37°C for 15 minutes. The reaction was terminated by the addition of 100 μl of BIOMOL® Green reagent and a subsequent incubation at room temperature for 30 minutes. Absorbance was measured using a spectrophotometer at 650 nm.</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

