

## Product Information

<b>Construct:</b>	PKCβ1 (GST-2-end)
<b>Concentration:</b>	0.10 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	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.
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	X06318
<b>MW:</b>	102 kDa
<b>Purity:</b>	80%
<b>Specific Activity:</b>	325 pmol/min/μg

Kinase activity was measured using a PKCtide peptide substrate (ERM<sub>2</sub>RPRKRQGSVRRRRV) diluted in distilled water at 1mg/ml.

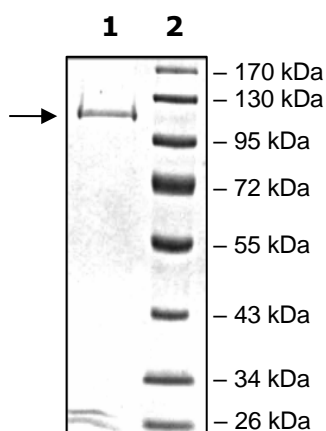
The protein kinase was diluted to 0.1 μg/μl in buffer containing 5 mM MOPS, pH 7.2, 2.5 mM β-glycero-phosphate, 5 mM MgCl<sub>2</sub>, 1 mM EGTA, 0.4 mM EDTA and 0.05 mM fresh DTT. Increasing amounts of the protein kinase were mixed with 0.38 mg/ml substrate and a sonicated PKC lipid activator in 20 μl final volume. The blank was determined from a "no substrate" sample. The reaction was initiated by addition of 5 μl of [33P]-ATP diluted in kinase buffer: 6ml kinase buffer containing 1 mCi [33P]-ATP, 0.25 mM ATP, 25 mM MOPS, pH 7.2, 12.5 mM β-glycero-phosphate, 25 mM MgCl<sub>2</sub>, 5 mM EGTA, 2 mM EDTA, and 0.25 mM fresh DTT.

After incubating for 30°C for 15 minutes, the reaction was 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

