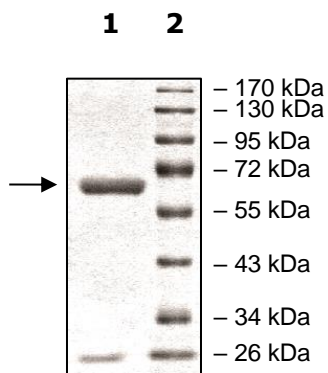


Product Information

Construct:	PKA α (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_002730
MW:	69 kDa
Purity:	≥90%
Specific Activity:	2,100 pmol/min/μg
Assay Conditions:	Kinase activity was measured using a CREBtide synthetic peptide substrate (KRREILSRPYSYR) at 1mg/ml. The protein kinase was diluted to 0.1 μg/μl in buffer containing 50ng/μl BSA and 5% glycerol. Increasing amounts of the protein kinase were mixed with 0.25 mg/ml substrate 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, 12.5 mM β-glycero-phosphate, 25 mM MgCl ₂ , 5 mM EGTA, 2 mM EDTA, 0.25 mM fresh DTT, pH 7.2. 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

