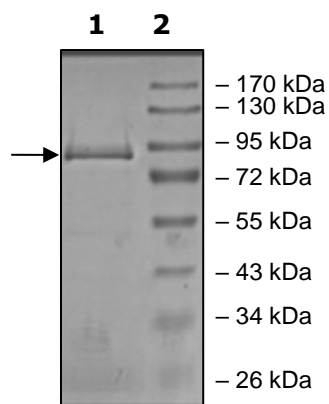


Product Information

Construct:	KAT5 (GST-Full length)
Concentration:	0.05 mg/ml
Species:	Human
Formulated In:	50 mM Tris-HCl, pH 7.5, 50 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_006388
MW:	88 kDa
Purity:	90%
Specific Activity:	30 pmol/min/μg
Assay Conditions:	Kinase activity was measured using Histone H4 Peptide substrate diluted in water to a final concentration of 1 mg/ml. Increasing amounts of the protein kinase were mixed with 0.25 mg/ml substrate to a 20 μl final volume. The reaction was initiated by addition of 5 μl of [Acetyl 3H]-CoA diluted in Acetyltransferase Assay Buffer, containing 50 mM Tris-HCL, pH 8.0, 0.1 mM EDTA, 5% glycerol and 2 mM fresh DTT for a final reaction volume of 25 μl. After incubating at 30°C for 15 minutes, the reaction was terminated by spotting 20 μl of the mixture onto phosphocellulose paper strips. Once fixed, strips were sequentially washed for at least three times in 50 mM Na ₂ HPO ₄ , pH 9.0 with gentle stirring. The blank was determined from a "no substrate" sample. 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

