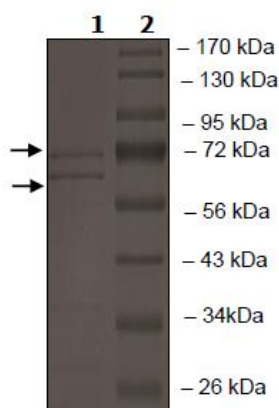


Product Information

Construct:	CDK14(GST-2-451(end))/Cyclin Y(GST-2-341(end))
Concentration:	0.10 mg/ml
Species:	Human
Formulated In:	50 mM Tris-HCl, pH 7.5, 50 mM NaCl, 10 mM glutathione, 0.25 mM DTT, 0.1 mM EDTA, 0.1 mM PMSF, 25% glycerol.
Expression System:	Co-expressed in Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	CDK14: NM_021395 Cyclin Y: NM_145012
MW:	CDK14: 72 kDa Cyclin Y: 64 kDa
Purity:	85%
Specific Activity:	≥2.5 pmol/min/μg
Assay Conditions:	5 mM MOPS, 2.5 mM β-glycerophosphate, 5 mM MgCl ₂ , 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 50 ng/ml BSA, 2 mM ATP, 200 μg of MBP (or Histone H1) as substrate. Add [33P]-ATP and incubate at 30°C for 15 minutes, then spot reaction on phosphocellulose paper, fix in 1% phosphoric.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

