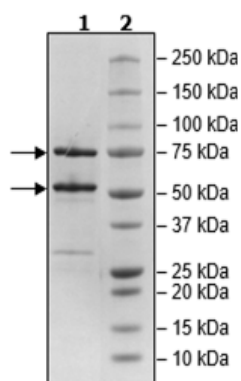


Product Information

Construct:	CDK1 (GST-Pre-2-297(end)) / CyclinA2 (GST-Pre-2-432(end))
Concentration:	0.84 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT
Expression System:	Coexpressed in Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	CDK1: NM_001786; CyclinA2: NM_001237
MW:	CDK1: 60 kDa; CyclinA2: 75 kDa
Purity:	≥90%
Specific Activity:	205 pmol/min/μg
Assay Conditions:	Assay was done according to the CDK1 Assay Kit (BPS Bioscience #79597) with various amounts of CDK1/CyclinA2.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

