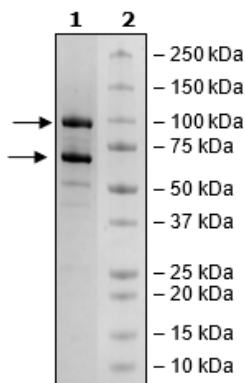


Product Information

Construct:	CDK12 (GST-696-1082) / CyclinK (GST-2-580(end))
Concentration:	0.33 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 3 mM DTT
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	CDK12: NM_016507; Cyclin K: NM_001099402
MW:	CDK12: 72 kDa; Cyclin K: 90 kDa
Purity:	≥90%
Specific Activity:	8 pmol/min/μg
Assay Conditions:	Assay was run according to CDK12/Cyclin K Kinase Assay Kit (BPS Bioscience #78298) in a Kinase buffer containing 2 mM DTT using 0.4 mg/ml CDK12/CyclinK substrate as a substrate with 10 μM ATP at 30°C for 90 min. The amount of ATP transferred from autophosphorylation as well as substrate phosphorylation was calculated using Kinase-Glo reagent (Promega).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CDK12/Cyclin K Activity

