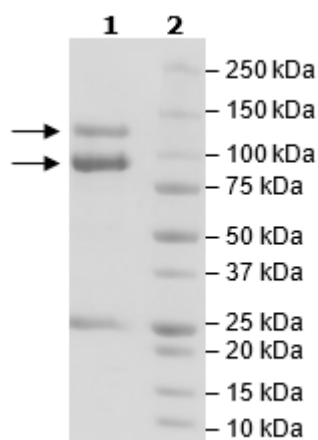


## Product Information

<b>Construct:</b>	CDK12 (GST-658-1490(end)) / CyclinK (GST-2-580(end))
<b>Concentration:</b>	0.36 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	CDK12: NM_016507; Cyclin K: NM_001099402
<b>MW:</b>	CDK12: 119 kDa; Cyclin K: 90 kDa
<b>Purity:</b>	80%
<b>Specific Activity:</b>	3 pmol/min/μg
<b>Assay Conditions:</b>	Assay was done according to CDK12/Cyclin K Kinase Assay Kit (BPS Bioscience #78298) with CDK12/Cyclin K titrated from 4000 ng/rxn to 0 ng/rxn.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

