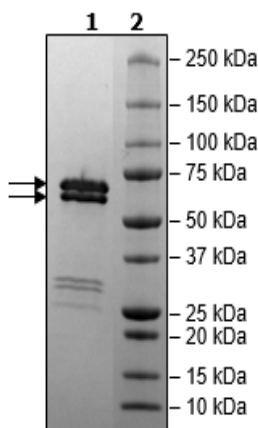


Product Information

<b>Construct:</b>	CDK16 (GST-107-496(end)) / CyclinY (GST-2-341(end))
<b>Concentration:</b>	0.93 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT and 20 mM glutathione
<b>Expression System:</b>	Coexpressed in 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>	CDK16: NM_006201; Cyclin Y: NM_145012
<b>MW:</b>	CDK16: 71 kDa; Cyclin Y: 66 kDa
<b>Purity:</b>	86%
<b>Specific Activity:</b>	≥35 pmol/min/μg
<b>Assay Conditions:</b>	Assay was done according to the CDK1 Assay Kit (BPS Bioscience #79597) with various amounts of CDK16.
<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

