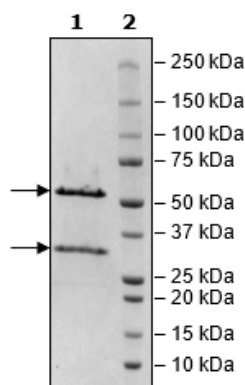


## Product Information

<b>Construct:</b>	CDK2 (1-298(end) / CyclinA2 (His-GST-174-432(end))
<b>Concentration:</b>	0.16 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	20 mM NaH <sub>2</sub> PO <sub>4</sub> /Na <sub>2</sub> HPO <sub>4</sub> , pH 8.0, 400 mM NaCl, 10% glycerol, 2 mM DTT, and 1 mM EDTA
<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>	CDK2: NM_001798; CyclinA2: NM_001237
<b>MW:</b>	CDK2: 34 kDa; CyclinA2: 57 kDa
<b>Purity:</b>	≥90%
<b>Specific Activity:</b>	203 pmol/min/μg
<b>Assay Conditions:</b>	Assay was done according to the CDK2 Assay Kit (BPS Bioscience #79599) with various amounts of CDK2/CyclinA2.
<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

