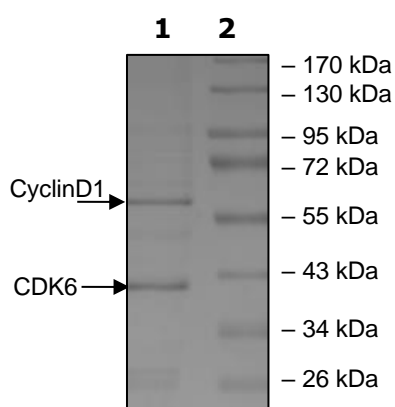


Product Information

Construct:	CDK6 (His-Full length) / CyclinD1 (GST-Full length)
Concentration:	0.10 mg/ml
Species:	Human
Formulated In:	50 mM Sodium Phosphate, pH 7.0, 300 mM NaCl, 150 mM imidazole, 0.1 mM PMSF, 0.2 mM DTT, 25% glycerol
Expression System:	Co-expressed Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	CDK6: NM_001259; CyclinD1: NM_053056
MW:	CDK6: 40 kDa; CyclinD1: 61 kDa
Purity:	80%
Specific Activity:	1.7 pmol/min/μg
Assay Conditions:	The activity of the complex was measured using ADP-Glo™ Kinase Assay kit (Promega; Cat# V9101) which quantifies the amount of ADP produced by phosphorylation. The ADP-Glo™ Reagent is added to terminate the reaction and quench the remaining ATP. The Kinase Detection Reagent is then added to convert ADP to ATP and to measure the newly converted ATP using a luciferase reaction. Assay: Kinase activity was measured using an Rb protein substrate. The reaction was initiated by mixing increasing amounts of the protein kinase with 25 μM ATP in 40 mM Tris-HCl, pH 7.4, 20 mM MgCl ₂ , 0.1 mg/ml BSA, 50 μM fresh DTT, and Rb substrate (0.04 mg/ml final). The reaction was terminated by addition of an equal volume of the ADP-Glo™ Reagent supplemented with 10mM MgCl ₂ , and the Kinase Detection Reagent was added. Phosphorylation was measured by detection of luminescence. The blank was determined from a “no kinase” sample.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

