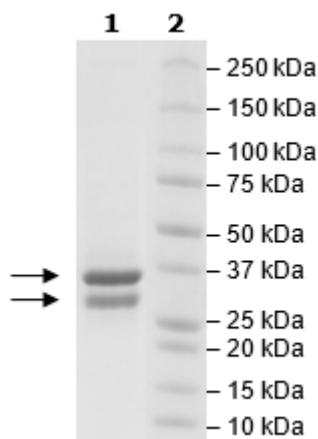


Product Information

Construct:	CDK6 (His-2-326(end)) / CyclinD3 (His-2-292(end))
Concentration:	1.76 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT and 20% glycerol
Expression System:	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; CyclinD3: NM_001760
MW:	CDK6: 38 kDa; CyclinD3: 33 kDa
Purity:	≥90%
Specific Activity:	50 pmol/min/μg
Assay Conditions:	Assay was done in a Kinase buffer using Histone H1 (0.1 mg/ml) as a substrate with 20 μM ATP at 30°C for 60 min. The amount of ATP transferred 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



CDK6/CyclinD3 Activity

