

CDK6/CyclinD3, Unactive

Full-length recombinant proteins expressed in Sf9 cells

Catalog # C35-14H

Lot # Q282-3

Product Description

Recombinant full-length human CDK6 and CyclinD3 were co-expressed by baculovirus in Sf9 insect cells using N-terminal His tags on both proteins. The gene accession numbers for CDK6 and CyclinD3 are [NM_001259](#) and [NM_001760](#), respectively.

Gene Aliases

PLSTIRE, MGC59692 /CCND3

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.2mM DTT, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

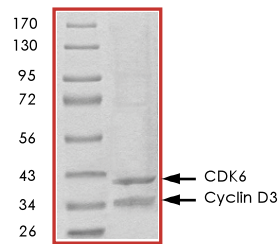
CDK6 is a member of the cyclin-dependent family of protein kinases that are important regulators of cell cycle progression. CDK6 activity is regulated by the D-type cyclins and members of the INK4 family of CDK inhibitors (1). The CDK6 kinase activity is detected in mid-G1 phase of the cell cycle and is responsible for the phosphorylation and regulation of the activity of tumor suppressor protein Rb. Although CDK6 and CDK4 can both phosphorylate multiple residues in the Rb protein, they do so with different residue selectivities in vitro; CDK6 phosphorylates Thr821 while CDK4 phosphorylates Thr826 on Rb protein (2).

References

- Meyerson, M. et al: Identification of G1 kinase activity for cdk6, a novel cyclin D partner. *Molec. Cell. Biol.* 14: 2077-2086, 1994.
- Takaki, T. et al: Preferences for phosphorylation sites in the retinoblastoma protein of D-type cyclin-dependent kinases, Cdk4 and Cdk6, in vitro. *J Biochem.* 2005 Mar; 137(3):381-6.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Purity



The purity of CDK6/CyclinD3 was determined to be **>90%** by densitometry, CDK6 approx. MW **40kDa** and CyclinD3 Approx. MW **35kDa**.

Upstream Active Kinases

CDK7/CyclinH1/MNAT1

Cat # C36-10H

CDK6/CyclinD3, Unactive

Full-length recombinant proteins expressed in Sf9 cells

Catalog Number C35-14H

Specific Lot Number Q282-3

Purity >90%

Concentration 0.2 $\mu\text{g}/\mu\text{l}$

Stability 1yr at -70°C from date of shipment

Storage & Shipping

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.