

## CyclinD1 Protein

Full length recombinant protein expressed in Sf9 cells

**Catalog # C81-30BG**

Lot # N044-2

### Product Description

Recombinant full-length human CyclinD1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_053056](#).

### Gene Aliases

BCL1, PRAD1, U21B31, D11S287E

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, and 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . 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

Cyclin D1 belongs to the highly conserved cyclin family, whose members are characterized by a dramatic periodicity in protein abundance throughout the cell cycle that functions as regulators of CDK kinases. Cyclin D1 is the regulatory subunit of a holoenzyme that phosphorylates and inactivates the RB protein and promotes progression through the G1-S phase of the cell cycle in a manner dependent on cyclin-dependent kinases, or CDKs. Cyclin D1 plays pivotal roles in the development of several human cancers, including parathyroid adenoma, breast cancer, colon cancer, lymphoma, melanoma, and prostate cancer (1). Cyclin D1 also coordinates mitochondrial function and nuclear DNA synthesis (2).

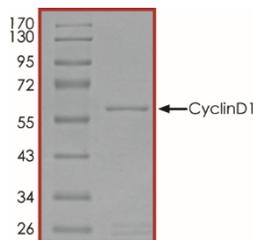
### References

1. Fu, M. et.al: Minireview: Cyclin D1: normal and abnormal functions. *Endocrinology* 145: 5439-5447, 2004.
2. Wang, C.et.al: Cyclin D1 repression of nuclear respiratory factor 1 integrates nuclear DNA synthesis and mitochondrial function. *Proc. Nat. Acad. Sci.* 103: 11567-11572, 2006.

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

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**

### Purity



The purity of CyclinD1 was determined to be **>90%** by densitometry.  
Approx. MW **61kDa**.

## CyclinD1 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog #	C81-30BG
Lot #	N044-2
Purity	>90%
Concentration	0.2 µg/µl
Stability	1yr at $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . 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.