

Cyclin T2 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # C86-30BG

Lot # K1623-4

Product Description

Full-length recombinant human Cyclin T2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_001241](#).

Gene Aliases

CCNT2; CYCT2

Formulation

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

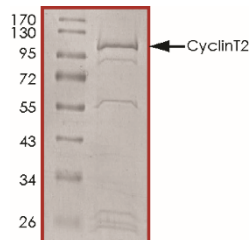
Scientific Background

Cyclin T2 belongs to the highly conserved cyclin family, whose members are characterized by a dramatic periodicity in protein abundance through the cell cycle which function as regulators of CDK kinases. Different cyclins exhibit distinct expression and degradation patterns which contribute to the temporal coordination of each mitotic event. Cyclin T2 and its kinase partner CDK9 were found to be subunits of the transcription elongation factor p-TEFb. The p-TEFb complex containing this cyclin was reported to interact with, and act as a negative regulator of human immunodeficiency virus type 1 (HIV-1) Tat protein(1). In mammalian cells, all 3 CDK9/cyclin T combinations strongly activated a CMV promoter(2).

References

1. Yang, Z.et.al: The 7SK small nuclear RNA inhibits the CDK9/cyclin T1 kinase to control transcription. Nature 414: 317-322, 2001.
2. Peng, J.et.al: Identification of multiple cyclin subunits of human P-TEFb. Genes Dev. 12: 755-762, 1998.

Purity



The purity of Cyclin T2 protein was determined to be **>70%** by densitometry. Approx. MW **102kDa**.

Cyclin T2 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	C86-30BG
Lot #	K1623-4
Purity	>70%
Concentration	0.1 µg/µl
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.

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

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