

CyclinA2 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # C78-30BG

Lot # N104-2

Product Description

Recombinant full-length human CyclinA2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_001237](#).

Gene Aliases

CCNA2; CCN1; CCNA

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM 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.

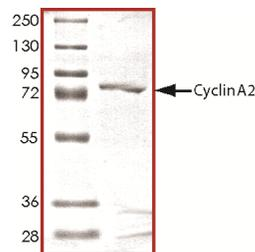
Scientific Background

CyclinA2 belongs to the highly conserved cyclin family whose members are characterized by a dramatic periodicity in protein abundance through the cell cycle and function as regulators of CDKs. CyclinA2 is an essential component of all embryonic and somatic cell cycles in mammals (1). CyclinA2 is highly expressed in the testis and certain myeloid leukemia cells (2). Cyclin A also plays a major role in the control of DNA replication (3). CyclinA2 binds and activates CDC2 or CDK2 and thus promotes both cell cycle G1/S and G2/M transitions.

References

1. Murphy, M.et.al: Delayed early embryonic lethality following disruption of the murine cyclin A2 gene. Nature Genet. 15: 83-86, 1997.
2. Yang, R.et.al: Characterization of a second human cyclin A that is highly expressed in testis and in several leukemic cell lines. Cancer Res. 57: 913-920, 1997.
3. Girard, F. et.al: Cyclin A is required for the onset of DNA replication in mammalian fibroblasts. Cell 67: 1169-1179, 1991.

Purity



The purity of CyclinA2 was determined to be **>75%** by densitometry. Approx. MW **78 kDa**.

CyclinA2 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	C78-30BG
Specific Lot Number	N104-2
Purity	>75%
Concentration	0.2 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.