

CDK11/CyclinC Protein

Full-length human recombinant protein expressed in Sf9 cells

Catalog # **C42-34G**

Lot # M154-1

Product Description

Recombinant full-length human CDK11/CyclinC was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession numbers for CDK11 and CyclinC are [NM_015076](#) and [NM_005190](#), respectively.

Gene Aliases

CDK11: CDC2L6; CDK8-like; bA346C16.3; KIAA1028

CyclinC: CCNC, CycC

Formulation

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

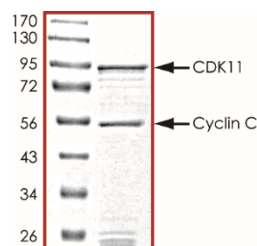
Scientific Background

CDK11/Cyclin C is a member of the p34Cdc2 protein kinase family which are known to be essential for eukaryotic cell cycle control. CDK11/Cyclin C gene complex may harbor 1 or more tumor suppressor genes affected by chromosome 1p36 modifications in neuroblastoma (1). CDK11/Cyclin C expression is regulated during murine embryogenesis and its activity is coordinately regulated with that of p34 (cdc2) during the cell cycle (2).

References

1. Lahti, J. M., et.al: Alterations in the PITSLRE protein kinase gene complex on chromosome 1p36 in childhood neuroblastoma. Nature Genet. 7: 370-375, 1994.
2. Bunnell, B.et.al: Elevated expression of a p58 protein kinase leads to changes in the CHO cell cycle. Proc. Nat. Acad. Sci. 87: 7467-7471, 1990. Note: Erratum: Proc. Nat. Acad. Sci. 88: 2612 only, 1991.

Purity



The purity of CDK11/CyclinC was determined to be **>90%** by densitometry, CDK11 approx. MW **92kDa**, and Cyclin H1 approx. MW **55kDa**.

CDK11/CyclinC Protein

Full-length human recombinant protein expressed in Sf9 cells

Catalog #	C42-34G
Lot #	M154-1
Purity	>90%
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.