

CDK11, Unactive

Recombinant full-length protein expressed in Sf9 cells

Catalog # C42-14G

Lot # S226-1

Product Description

Recombinant full-length human CDK11 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The CDK11 gene accession number is [NM_015076](#).

Gene Aliases

CDK19; CDC2L6; CDK8-like; bA346C16.3; KIAA1028

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.

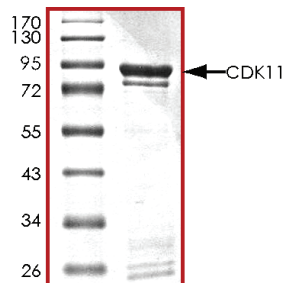
Scientific Background

CDK11 is a member of the cyclin-dependent kinase 11 family which is known to be essential for eukaryotic cell cycle control and it is in close proximity to CDC2L2, a nearly identical gene in the same chromosomal region. CDK11 is a negative regulator of normal cell cycle progression and its expression is regulated during murine embryogenesis, and its activity is coordinately regulated with that of p34 (cdc2) during the cell cycle (1). CDK11 gene may harbor 1 or more tumor suppressor genes that affected by chromosome 1p36 modifications in neuroblastoma (2). CDK11 could be cleaved by caspases and was demonstrated to play roles in cell apoptosis.

References

1. 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.
2. 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.

Purity



The purity of CDK11 was determined to be **>90%** by densitometry, approx. MW **92 kDa**.

CDK11, Unactive

Recombinant full-length human protein expressed in Sf9 cells

Catalog Number	C42-14G
Specific Lot Number	S226-1
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.