

## CyclinK Protein

Full length recombinant protein expressed in Sf9 cells

**Catalog # C84-30G**

Lot # N092-2

### Product Description

Recombinant full-length human CyclinK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC015935](#).

### Gene Aliases

CCNK, CPR4, MGC9113

### 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.

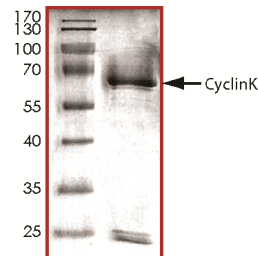
### Scientific Background

CyclinK is a member of the cyclin family which can regulate transcription through their association with and activation of cyclin-dependent kinases (CDK). The CDK/CyclinK complex can phosphorylate the C-terminal domain (CTD) of the large subunit of RNA polymerase II (1). CyclinK is associated with potent CTD kinase and CDK kinase activities in vitro and coimmunoprecipitates with POLR2A. CyclinK is a new member of the transcription cyclin family which may play a dual role in regulating CDK and RNA polymerase II activities (2).

### References

1. Edwards, M. C. et.al: Human cyclin K, a novel RNA polymerase II-associated cyclin possessing both carboxy-terminal domain kinase and Cdk-activating kinase activity. *Molec. Cell. Biol.* 18: 4291-4300, 1998.
2. Lin X: P-TEFb containing cyclin K and Cdk9 can activate transcription via RNA. *J Biol Chem.* 2002 May 10;277(19):16873-8.

### Purity



The purity of CyclinK was determined to be **>85%** by densitometry. Approx. MW **67 kDa**.

## CyclinK Protein

Full length recombinant protein expressed in Sf9 cells

Catalog #	C84-30G
Lot #	N092-2
Purity	>85%
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](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.**