

## CDK13/CyclinK Protein

Recombinant human protein expressed in Sf9 cells

**Catalog # C25-35BG**

Lot # G1404-1

### Product Description

Recombinant human CDK13 (516-end) /Cyclin K complex was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins. The gene accession number for CDK13 and CyclinK are [NM\\_003718](#) and [BC015935](#), respectively.

### Gene Aliases

CDK13: CDC2L; CDC2L5; CHED; hCDK13  
CyclinK: 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.

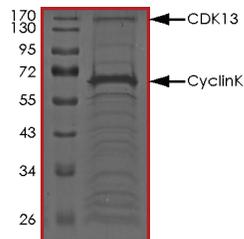
### Scientific Background

CDK13 or cyclin-dependent kinase 13 is a member of the cyclin-dependent serine/threonine protein kinase family which are well known for their essential roles as master switches in cell cycle control. CDK13 forms a complex with cyclin K and is predicted to play a role in gene regulation (1). CDK13 may play a role in mRNA processing and is thought to be involved in regulation of hematopoiesis (2). The CYCK-CDK12 and CYCK-CDK13 complexes display unique function in cell cycle and gene regulation.

### References

- Blazek, D.et.al: The cyclin K/Cdk12 complex maintains genomic stability via regulation of expression of DNA damage response genes. *Genes Dev.* 25: 2158-2172, 2011.
- Lapidot-Lifson, Y.et.al: Cloning and antisense oligodeoxynucleotide inhibition of a human homolog of cdc2 required in hematopoiesis. *Proc. Nat. Acad. Sci.* 89: 579-583, 1992.

### Purity



The purity of CDK13/CyclinK was determined to be **>70%** by densitometry, CDK13 approx. MW **160kDa** and CyclinK Approx. MW **67kDa**.

## CDK13/CyclinK Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	C25-35BG
Specific Lot #	G1404-1
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](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.**