

## CyclinH Protein

Full length recombinant protein expressed in insect cells

**Catalog # C83-30H**

Lot # O797-3

### Product Description

Recombinant full length human CyclinH was expressed in insect cells using an N-terminal His tag. The gene accession number is [NM\\_001239](#).

### Gene Aliases

CCNH, CAK, p34, p37

### Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 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.

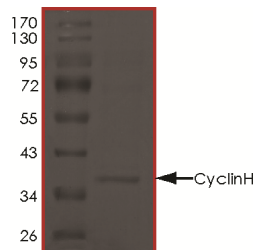
### Scientific Background

Cyclin H is related to the highly conserved cyclin family. Cyclin family members are characterized by a dramatic periodicity in protein abundance through the cell cycle. Cyclins function as regulators of CDK kinases. Cyclin H forms a complex with CDK7 kinase and the ring finger protein MAT1 which is able to phosphorylate CDK2 and CDC2 kinases and thus act as a CDK-activating kinase (CAK) (1). Cdk-activating kinase complex is a component of human transcription factor TFIH and plays an important role in the processes of transcription, DNA repair, and cell cycle progression (2).

### References

1. Eki, T. et.al: Mapping of the human genes encoding cyclin H (CCNH) and the CDK-activating kinase (CAK) assembly factor MAT1 (MNAT1) to chromosome bands 5q13.3-q14 and 14q23, respectively. Genomics 47: 115-120, 1998.
2. Shiekhhattar, R. et.al: Cdk-activating kinase complex is a component of human transcription factor TFIH. Nature 374: 283-287, 1995.

### Purity



The purity of CyclinH was determined to be **>90%** by densitometry. Approx. MW **39kDa**.

## CyclinH Protein

Full length recombinant protein expressed in insect cells

Catalog Number C83-30H

Specific Lot Number O797-3

Purity	>90%
Concentration	0.1µg/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.**