

Catalog #	Aliquot Size
C38-34G-20	20 µg
C38-34G-50	50 µg

CDK8/CyclinC Protein

Full-length human recombinant protein expressed in Sf9 cells

Catalog # C38-34G

Lot # M082-1

Product Description

Recombinant full-length human CDK8/CyclinC was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession numbers for CDK8 and CyclinC are [BC069634](#) and [NM_005190](#), respectively.

Gene Aliases

CDK8: K35, MGC126074, MGC126075

CyclinC: CCNC, CycC

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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

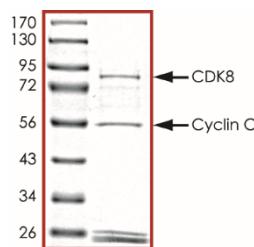
CDK8 is a member of the CDK family which are important regulators of cell cycle progression. CDK8 associates with Cyclin C and is a component of the RNA polymerase II holoenzyme complex that phosphorylates the carboxy-terminal domain (CTD) of the largest subunit of RNA polymerase II (1). CDK8 regulate transcription by targeting the CDK7/cyclin H subunits of the general transcription initiation factor IIH (TFIIH). Inhibition of CDK8 expression inhibits proliferation in colon cancer cells characterized by high levels of CDK8 and beta-catenin hyperactivity. CDK8 kinase activity is necessary for beta-catenin-driven transformation and for expression of several beta-catenin transcriptional targets (2).

References

1. Akoulitchev S, et al: TFIIH is negatively regulated by cdk8-containing mediator complexes. *Nature* 407: 102-106, 2000.
2. Firestein R, et al: CDK8 is a colorectal cancer oncogene that regulates beta-catenin activity. *Nature* 455: 547-551, 2008.

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

Purity



The purity of CDK8/CyclinC was determined to be >70% by densitometry, CDK8 approx. MW **84kDa**, and Cyclin H1 approx. MW **55kDa**.

CDK8/CyclinC Protein

Full-length human recombinant protein expressed in Sf9 cells

Catalog #	C38-34G
Lot #	M082-1
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.

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.