

CDK7, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog # C36-14G

Lot # S334-1

Product Description

Recombinant full-length human CDK7 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tags. The gene accession number is [NM_001799](#).

Gene Aliases

CDK7: CAK1, STK1, CDKN7, p39MO15

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.

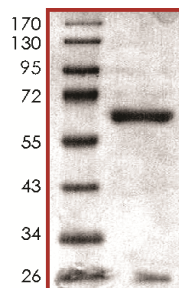
Scientific Background

CDK7 gene is a member of the cyclin-dependent protein kinase family that is important regulators of cell cycle progression (1). CDK7 forms a trimeric complex with cyclin H and MAT1, which functions as a CDK-activating kinase (CAK). CDK7 is an essential component of the transcription factor TFIIH that is involved in transcription initiation and DNA repair. CDK7 is thought to serve as a direct link between the regulation of transcription and the cell cycle (2).

References

1. Fisher, R. P.; A novel cyclin associates with MO15/CDK7 to form the CDK-activating kinase. *Cell* 78: 713-724, 1994.
2. Larochelle, S. Requirements for Cdk7 in the assembly of Cdk1/cyclin B and activation of Cdk2 revealed by chemical genetics in human cells. *Molec. Cell* 25: 839-850, 2007.

Purity



The purity of CDK7 was determined to be **>95%** by densitometry. Approx. MW **66 kDa**.

CDK7, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	C36-14G
Specific Lot Number	S334-1
Purity	>95%
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.