

CDKN3 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog # C01-34G

Lot # K125-1

Product Description

Recombinant full-length human CDKN3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_005192](#).

Gene Aliases

CDI1, CIP2, FLJ25787, KAP, KAP1, MGC70625

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

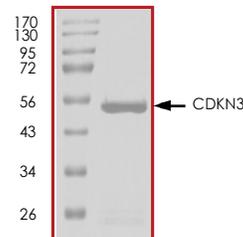
CDKN3 is a member of the dual specificity protein phosphatase family that acts as a cyclin-dependent kinase inhibitor (1). CDKN3 has been shown to interact with and dephosphorylate specifically the CDK2 kinase thereby preventing its activation. CDKN3 has been reported to be deleted, mutated, or overexpressed in several kinds of cancers. Increased expression of CDKN3 leads to increased levels of kinase-associated phosphatase activity that inhibits the G(1)/S transition of the cell cycle by dephosphorylating the cyclin-dependent kinases (2).

References

1. Niculescu, M.D. et al : Choline availability modulates human neuroblastoma cell proliferation and alters the methylation of the promoter region of the cyclin-dependent kinase inhibitor 3 gene. *J Neurochem.* 2004, 89(5):1252-9.
2. Yeh, C.T. et al: Abolishment of the interaction between cyclin-dependent kinase 2 and Cdk-associated protein phosphatase by a truncated KAP mutant. *Biochem. Biophys. Res. Commun.* 2003; 305(2):311-4

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 CDKN3 was determined to be **>90%** by densitometry. Approx. MW **51kDa**.

CDKN3 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	C01-34G
Specific Lot Number	K125-1
Purity	>90%
Concentration	0.2 µg/µl
Stability	1 yr 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.