

CyclinO Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # C85-30G

Lot # O869-1

Product Description

Recombinant full-length human CyclinO was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_021147](#).

Gene Aliases

CCNO; CCNU; UDG2

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.

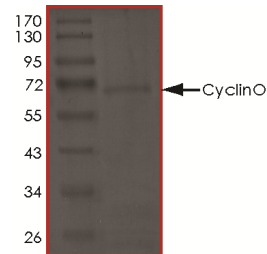
Scientific Background

Cyclin O belongs to the highly conserved cyclin family. These family members are characterized by a dramatic periodicity in protein abundance through the cell cycle. Cyclin O plays an important role in the processes of transcription. Cyclin O also plays an important role in DNA repair and cell cycle progression. Cyclin O localizes exclusively in the nucleus (1). Cyclin O shows significant uracil-DNA glycosylase activity in which the activity of in vitro translated UDG2 was not as robust as that of the mitochondrial isoform of UNG (2).

References

1. Caradonna, S. et.al: Affinity purification and comparative analysis of two distinct human uracil-DNA glycosylases. *Exp. Cell Res.* 222: 345-359, 1996.
2. Muller, S. J. et.al: Cell cycle regulation of a human cyclin-like gene encoding uracil-DNA glycosylase. *J. Biol. Chem.* 268: 1310-1319, 1993.

Purity



The purity of CyclinO was determined to be **>95%** by densitometry. Approx. MW **69kDa**.

CyclinO Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	C85-30G
Specific Lot Number	O869-1
Purity	>90%
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
www.signalchem.com

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