

UHK1 (KIS) Protein

Recombinant full-length protein expressed in Sf9 cells

Catalog # K05-30G

Lot # X681-4

Product Description

Recombinant full-length human UHK1 (KIS) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_144624](#).

Gene Aliases

UHK1; KIS; KIST

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.

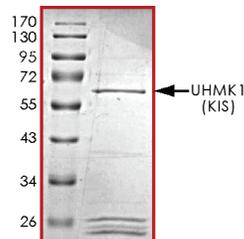
Scientific Background

UHK1(KIS) or U2AF homology motif (UHM) kinase 1 is a serine/threonine protein kinase that promotes cell cycle progression through G1 phase. UHK1 phosphorylates the cyclin-dependent kinase inhibitor 1B (p27Kip1) and this results in its export to the nucleus where it undergoes degradation. UHK1 is a growth factor-dependent nuclear kinase which can regulate cell cycle progression (1). Elevated levels of UHK1 protein in leukemia cells has been shown to promote cell cycle progression. UHK1 has also been shown to function in the adult nervous system where it is highly expressed and the gene has been associated with schizophrenia (2).

References

1. Boehm. et.al: A growth factor-dependent nuclear kinase phosphorylates p27(Kip1) and regulates cell cycle progression. EMBO J. 21: 3390-3401, 2002.
2. Bieche. et.al: Quantitative RT-PCR reveals a ubiquitous but preferentially neural expression of the KIS gene in rat and human. Molec. Brain Res. 114: 55-64, 2003.

Purity



The purity of UHK1 (KIS) was determined to be **>70%** by densitometry, approx. MW **65 kDa**.

UHK1 (KIS) Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog	K05-30G
Lot #	X681-4
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.

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.