

HEMK1 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog # H311-30G

Lot # L1882-4

Product Description

Recombinant full-length human HEMK1 protein was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The HEMK1 gene accession number is [NM_016173](#).

Gene Aliases

HEMK; MTQ1

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.

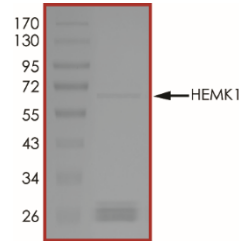
Scientific Background

HEMK1 belongs to the HemK methyltransferase family, members of which are found in genomes from bacteria to humans (1). HEMK1 methylates the glutamine residue in the GGQ motif of HMRF1L, a human mitochondrial release factor, and plays an important regulatory role in translation termination process (2).

References

1. Bujnicki, J.M., et al. Is the HemK family of putative S-adenosylmethionine-dependent methyltransferases a "missing" zeta subfamily of adenine methyltransferases? A hypothesis. IUBMB Life. 48(3):247-9, 1999.
2. shizawa, T., et al. The human mitochondrial translation release factor HMRF1L is methylated in the GGQ motif by the methyltransferase HMPmC. Biochem Biophys Res Commun. 373(1):99-103, 2008.

Purity



The purity of HEMK1 was determined to be >70% by densitometry, approx. MW **65 kDa**.

HEMK1 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog #	H311-30G
Lot #	L1882-4
Purity	>70%
Concentration	0.05 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.

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.