

KDM4C Protein

Recombinant protein expressed in Sf9 cells

Catalog # K424-31CG

Lot # J606-1

Product Description

Recombinant human KDM4C (1-460) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC104861](#).

Gene Aliases

KDM4C, GASC1, JHDM3C, JMJD2C

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.

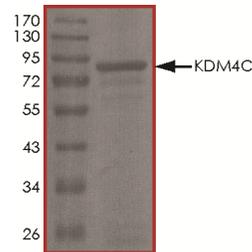
Scientific Background

KDM4C or lysine-specific demethylase 4C functions as a trimethylation-specific demethylase, converting specific trimethylated histone residues to the dimethylated form (1). KDM4C is a member of the Jumonji domain 2 (JMJD2) family and contains one JmjC domain, one JmjN domain, two PHD-type zinc fingers, and two Tudor domains. KDM4C may contribute to tumor development and inhibition of KDM4C expression leads to decreased cell proliferation. KDM4C has been shown to play an important role in the development and/or progression of various types of cancer, including esophageal squamous cell carcinoma(2).

References

1. Cloos, P. A. C.et.al: The putative oncogene GASC1 demethylates tri- and dimethylated lysine 9 on histone H3. *Nature* 442: 307-311, 2006.
2. Yang, Z.-Q.et.a: Identification of a novel gene, GASC1, within an amplicon at 9p23-24 frequently detected in esophageal cancer cell lines. *Cancer Res.* 60: 4735-4739, 2000.

Purity



The purity of KDM4C was determined to be **>95%** by densitometry. Approx. MW **82 kDa**.

KDM4C Protein

Recombinant protein expressed in Sf9 cells

Catalog Number	K424-31CG
Specific Lot Number	J606-1
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.

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.