

KDM4A Protein

Recombinant human protein expressed in Sf9 cells

Catalog # K424-31G

Lot # 1291-2

Product Description

Recombinant human KDM4A (1-886) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC002558](#).

Gene Aliases

JMJD2A; JHDM3A; KIAA0677

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.

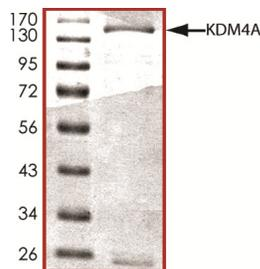
Scientific Background

KDM4A or lysine (K)-specific demethylase 4A is a member of the Jumonji domain 2 (JMJD2) family that encodes a protein containing a JmjN domain, a JmjC domain, a JD2H domain, two TUDOR domains, and two PHD-type zinc fingers. KDM4A functions in euchromatin to remove histone methylation marks that are associated with active transcription (1). KDM4A also act as a trimethylation-specific demethylase those converting specific trimethylated histone residues to the dimethylated form, and as a transcriptional repressor (2).

References

1. Klose, R. J. et.al: The transcriptional repressor JHDM3A demethylates trimethyl histone H3 lysine 9 and lysine 36. *Nature* 442: 312-316, 2006.
2. Whetstine, J. R. et.al: Reversal of histone lysine trimethylation by the JMJD2 family of histone demethylases. *Cell* 125: 467-481, 2006.

Purity



The purity of KDM4A was determined to be **>90%** by densitometry. Approx. MW **150kDa**.

KDM4A Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number	K424-31G
Specific Lot Number	1291-2
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.