

KDM1B Protein

Full length mouse recombinant protein expressed in Sf9 cells

Catalog # K421-30BG

Lot # P1749-3

Product Description

Recombinant full-length mouse KDM1B was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_172262](#).

Gene Aliases

4632428N09Rik; AI482520; Aof1

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.

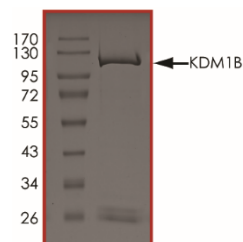
Scientific Background

KDM1B encodes Lysine-specific histone demethylase 1B that demethylates 'Lys-4' of histone H3, a specific tag for epigenetic transcriptional activation, thereby acting as a corepressor. It specifically associates with the coding region of its target genes. Removal of endogenous KDM1B promotes an increase in H3K4me2 levels and concurrent decrease in H3K9me2 levels, with a consequent down-regulation of targeted gene transcription.

References

1. Fang R, et al. Human LSD2/KDM1b/AOF1 regulates gene transcription by modulating intragenic H3K4me2 methylation. Mol Cell. 2010 Jul 30;39(2):222-33.
2. <http://www.uniprot.org/uniprot/Q8NB78>.

Purity



The purity of KDM1B was determined to be **>85%** by densitometry. Approx. MW **~118kDa**.

KDM1B Protein

Full length mouse recombinant protein expressed in Sf9 cells

Catalog Number

K421-30BG

Lot #

P1749-3

Purity

>85%

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.