

KDM5B Protein

Recombinant protein expressed in Sf9 cells

Catalog # K425-31BG

Lot # F600-3

Product Description

Recombinant mouse KDM5B (1-753) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC057318](#).

Gene Aliases

Jarid1b;PUT1;Plu1;PLU-1;Rb-Bp2; AW556288; RBBP2H1A; mKIAA4034

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.

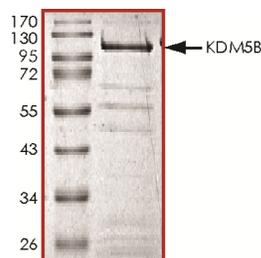
Scientific Background

KDM5B is a lysine (K)-specific demethylase 5B which is highly expressed in testis and shows the lowest expression in skeletal muscle (1). KDM5B is a histone H3 lysine 4 (H3K4) demethylase up-regulated in prostate cancer and associated with androgen receptor where it regulated its transcriptional activity in reporter gene assays (2). Overexpression of KDM5B in HeLa cells results in loss of tri-, di-, and monomethyl H3K4, but does not alter the methylation status of H3K9 or H3K36. A KDM5B containing highly conserved putative DNA/chromatin binding motif is specifically up-regulated in breast cancer.

References

1. Kashuba, V. et.al: Isolation and chromosomal localization of a new human retinoblastoma binding protein 2 homologue 1a (RBBP2H1A). *Europ. J. Hum. Genet.* 8: 407-413, 2000.
2. Xiang, Y. et.al: JARID1B is a histone H3 lysine 4 demethylase up-regulated in prostate cancer. *Proc. Nat. Acad. Sci.* 104: 19226-19231, 2007.

Purity



The purity of KDM5B was determined to be **>90%** by densitometry. Approx. MW **112kDa**.

KDM5B Protein

Recombinant protein expressed in Sf9 cells

Catalog Number	K425-31BG
Specific Lot Number	F600-3
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.