

KDM4B Protein

Recombinant protein expressed in Sf9 cells

Catalog # K424-31BG

Lot # F566-1

Product Description

Recombinant mouse KDM4B (1-725) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_172132.2](#).

Gene Aliases

Jmjd2b; mKIAA0876; 4732474L06Rik

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

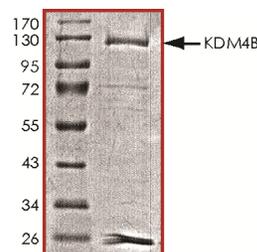
KDM4B or lysine (K)-specific demethylase 4B, which is also known as JMJD2B, contains a JmjN domain, a JmjC domain, a JD2H domain, 2 TUDOR domains, and a bipartite nuclear localization signal that overlaps the C-terminal part of the second TUDOR domain. KDM4B plays an essential role in human carcinogenesis through positive regulation of cyclin-dependent kinase 6 (1). KDM4B functions as a co-factor of estrogen receptor in breast cancer proliferation and mammary gland development (2). KDM4B is regulated by both ER α and HIF-1 α , drives breast cancer cell proliferation in normoxia and hypoxia, and epigenetically regulates the expression of cell cycle genes such as CCND1, CCNA1, and WEE1.

References

1. Toyokawa G. et.al: The histone demethylase JMJD2B plays an essential role in human carcinogenesis through positive regulation of cyclin-dependent kinase 6. 2011 Dec.
2. Kawazu M. et.al: Histone demethylase JMJD2B functions as a co-factor of estrogen receptor in breast cancer proliferation and mammary gland development, 2011 Mar 18.

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

Purity



The purity of KDM4B was determined to be **>75%** by densitometry. Approx. MW **125kDa**.

KDM4B Protein

Recombinant protein expressed in Sf9 cells

Catalog Number	K424-31BG
Specific Lot Number	F566-1
Purity	>75%
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.

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.