

Anti-KDM4B

Rabbit Polyclonal Antibody

Catalog # K424-63BR

Lot # O2121-24

Cited Applications

WB, ELISA

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the KDM4B protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

KDM4B antibody was raised against a 16 amino acid synthetic peptide from near the center of human KDM4B

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

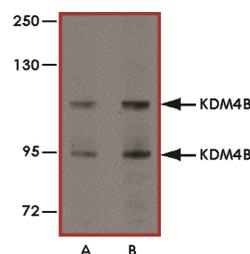
Scientific Background

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.

Sample Data



Western blot analysis of KDM4B in A20 cell lysate with KDM4B antibody at (A) 1 and (B) 2 µg/ml.

Anti-KDM4B

Rabbit Polyclonal Antibody

Catalog Number

K424-63BR

Specific Lot Number

O2121-24

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

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.