

Anti-KDM6B

Rabbit Polyclonal Antibody

Catalog # K426-63BR

Lot # O2361-31

Cited Applications

ELISA, WB, ICC, IF

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

Specificity

Recognizes the KDM6B protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

KDM6B antibody was raised against a 16 amino acid synthetic peptide from near the amino terminus of human KDM6B

Formulation

PBS + 0.02% sodium azide.

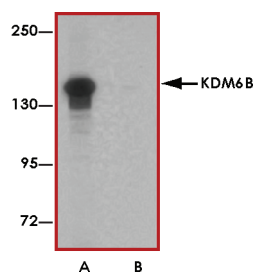
Stability

1yr at -20°C from date of shipment.

References

1. Lai F et al: cDNA cloning and genomic structure of three genes localized to human chromosome band 5q31 encoding potential nuclear proteins. *Genomics*2000; 70:123-30.
2. Agger K et al: UTX and JMJD3 are histone H3K27 demethylases involved in HOX gene regulation and development. *Nature* 2007; 449:731-4.
3. De Santa F et al: The histone H3 lysine-27 demethylase Jmjd3 links inflammation to inhibition of polycomb-mediated silencing. *Cell* 2007; 130:1083-94
4. Sola S et al: p53 interaction with JMJD3 results in its nuclear distribution during mouse neural stem cell differentiation. *PLoS One* 2011; 6:e18421.

Sample Data



Western blot analysis of KDM6B in K562 cell lysate with KDM6B antibody at 0.5 µg/mL in (A) the absence and (B) the presence of blocking peptide.



Immunocytochemistry of KDM6B in K562 cells with KDM6B antibody at 2.5 µg/mL.

Anti-KDM6B

Rabbit Polyclonal Antibody

Catalog Number

K426-63BR

Specific Lot Number

O2361-31

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.