

Anti-KDM1A

Rabbit Polyclonal Antibody

Catalog # K421-63R

Lot # O2109-49

Cited Applications

WB, ELISA

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

Specificity

Recognizes the KDM1A protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

KDM1A antibody was raised against a 17 amino acid synthetic peptide from near the amino terminus of human KDM1A

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

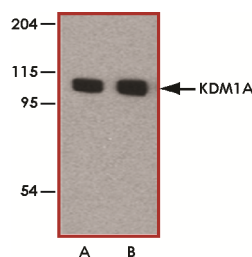
Scientific Background

KDM1A or lysine (K)-specific demethylase 1A is a nuclear protein containing a SWIRM domain, a FAD-binding motif, and an amine oxidase domain that is a component of several histone deacetylase complexes and act as a histone demethylase. (1) H3K4 histone demethylase activity of KDM1A is partly responsible for the repressive activity of TAL1 and restricts TAL1 function in hematopoiesis (1). KDM1A plays an essential role for CoREST in demethylation of H3K4 both in vitro and in vivo (2).

References

1. Hu, X. et.al : LSD1-mediated epigenetic modification is required for TAL1 function and hematopoiesis. Proc. Nat. Acad. Sci. 106: 10141-10146, 2009.
2. Lee, M. G. et.al : An essential role for CoREST in nucleosomal histone 3 lysine 4 demethylation. Nature 437: 432-435, 200.

Sample Data



Western blot analysis of KDM1A in P815 cell lysate with KDM1A antibody at (A) 1 and (B) 2 µg/ml.

Anti-KDM1A

Rabbit Polyclonal Antibody

Catalog Number

K421-63R

Specific Lot Number

O2109-49

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.