

KDM1B polyclonal antibody

Other names: AOF1, LSD2, C6orf193

Cat. No. C15310248 (CS-PA028-100)

Type: Polyclonal

Source: Rabbit

Lot #: A1224-004

Size: 100 µl

Concentration: not determined

Specificity: Human: positive / Other species: not tested

Purity: Whole antiserum from rabbit containing 0.05% azide.

Storage: Store at -20°C; for long storage, store at -80°C.
Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Description: Polyclonal antibody raised in rabbit against human KDM1B (lysine (K)-specific demethylase 1B) using two KLH-conjugated synthetic peptides containing a sequence from the N-terminal and the central region of the protein, respectively.

Applications

	Suggested dilution	Results
ELISA	1:500	Fig 1

*The optimal dilution for other applications should be determined by the end user. For WB we suggest starting with a 1:1,000 dilution

Target description

KDM1B (UniProtKB/Swiss-Prot entry Q8NB78) is a histone demethylase that demethylates lysine 4 of histone H3, a specific tag for epigenetic transcriptional activation, thereby acting as a corepressor. KDM1B is able to demethylate both mono- and di-methylated lysine 4 of histone H3, but has no effect on tri-methylated lysine 4, mono-, di- or tri-methylated lysine 9, mono-, di- or tri-methylated lysine 27, mono-, di- or tri-methylated lysine 36 of histone H3, or on mono-, di- or tri-methylated lysine 20 of histone H4. KDM1B is also required for de novo DNA methylation of a subset of imprinted genes during oogenesis. It acts by oxidizing the substrate by FAD to generate the corresponding imine that is subsequently hydrolyzed.

Results

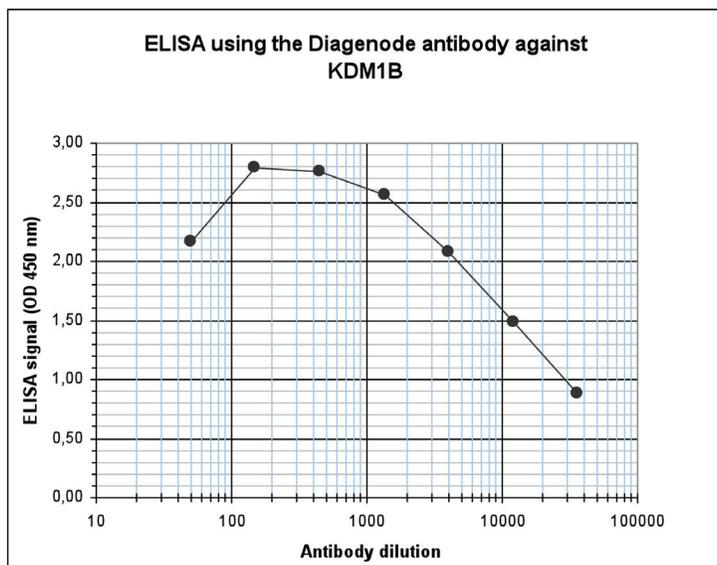


Figure 1. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human KDM1B (cat. No. CS-PA028-100). The plates were coated with the peptides used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:14,250.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: September 5, 2012