

PRODUCT NAME		
LSD1 polyclonal antibody		
Other names: BHC110, AOF2, EC1, KDM1		
Cat. No. C15410125 (pAb-125-050)	Type: Polyclonal	Size: 50 µg /75 µl
Lot #: A56-001	Source: Rabbit	Concentration: 0.67 µg/µl

Description: Polyclonal antibody raised in rabbit against human LSD1 (Lysine-specific demethylase 1), using a KLH-conjugated synthetic peptide containing a sequence from the N-terminal part of the protein.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:100	Fig 1
Western blotting	1:1,000	Fig 2, Fig 3
Immunoprecipitation	2 µg/IP	Fig 3

Purity: Affinity purified polyclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300.

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.

Last data sheet update: March 4, 2010

Target description

LSD1 (lysine specific demethylase 1, UniProt/Swiss-Prot entry O60341) is a component of the histone demethylase complex which uses FAD as a prosthetic group. LSD1 may have a dual effect on gene transcription. As it demethylates the mono- and dimethylated 'Lys-4' of histone H3, which are associated with transcriptional activation, LSD1 can act as a repressor of gene expression. However, LSD1 is also capable of demethylating 'Lys-9' of histone H3, a specific tag for epigenetic transcriptional repression, thereby leading to activation of androgen receptor target genes. As a consequence, LSD1 is involved in different processes such as embryonic development, cell differentiation and proliferation, and stem and cancer cell biology.

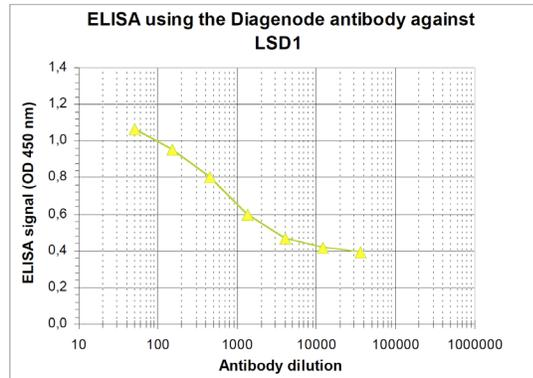


Figure 1
Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human LSD1 (Cat. No. pAb-125-050). The plates were coated with the peptide 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 1:2,500.



Figure 2
Western blot analysis using the Diagenode antibody directed against LSD1

Western blot was performed using respectively 150 µg (lane 1) or 50 µg (lane 2) nuclear extracts from HeLa cells and the Diagenode antibody against LSD1 (Cat. No. pAb-125-050) diluted 1:1,000. The location of the protein of interest is indicated on the left.

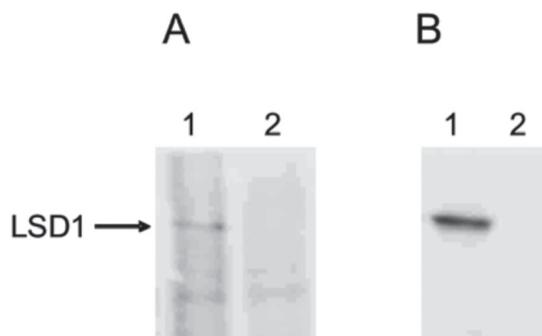


Figure 3

Immunoprecipitation and Western blot analysis using the Diagenode antibody directed against LSD1

293T cells were transiently transfected with an expression vector for HA-tagged LSD1 (lanes 1) or with an empty vector (lanes 2).

Figure 3A: Whole cell extracts were analysed by Western blot using the Diagenode antibody against human LSD1 (Cat. No. pAb-125-050), diluted 1:1,000.

Figure 3B: Whole cell extracts were immunoprecipitated with 2 µg of the Diagenode antibody against LSD1. Immunoprecipitates were analysed by Western blot as described above.