

LSD1 polyclonal antibody

Other names: BHC110, AOF2, EC1, KDM1

Cat. No. C15310028

Type: Polyclonal

Source: Rabbit

Lot #: 001

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 LSD1 (Lysine-specific demethylase 1), using the full length recombinant protein.

Applications

	Suggested dilution	Results
Western blotting	1:1,000	Fig 1, 2

Target description

LSD1 (lysine specific demethylase 1, UniProt/Swiss-Prot entry O60341) is a component of the histone demethylase complex that 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. LSD1 therefore mediates different processes such as embryonic development, cell differentiation and proliferation, stem and cancer cell biology.

Results

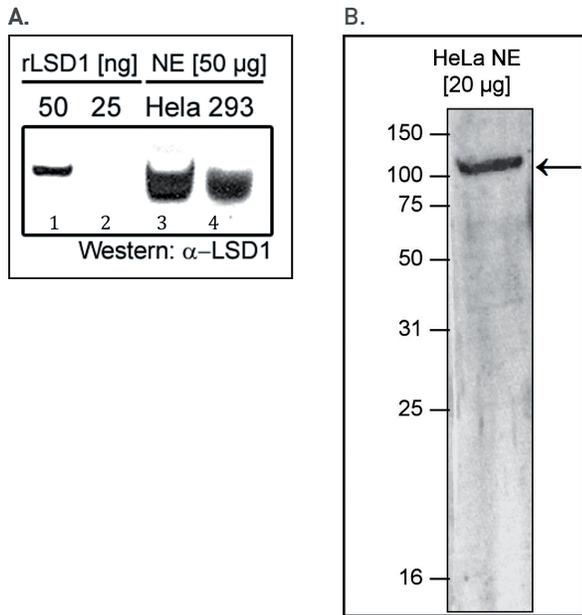


Figure 1. Western blot analysis using the Diagenode antibody directed against LSD1

Figure 1A: Western blot was performed with the Diagenode antibody against LSD1 (Cat. No. C15310028) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk on purified recombinant LSD1 protein (50 and 25 ng, lane 1 and 2) and on nuclear extracts (50 µg) from HeLa and HEK293 cells (lane 3 and 4, respectively).

Figure 1B: Western blot was performed as described above using nuclear extracts from HeLa cells (HeLa NE, 20 µg). The molecular weight marker (in kDa) is shown on the left; the location of the protein of interest is indicated on the right.

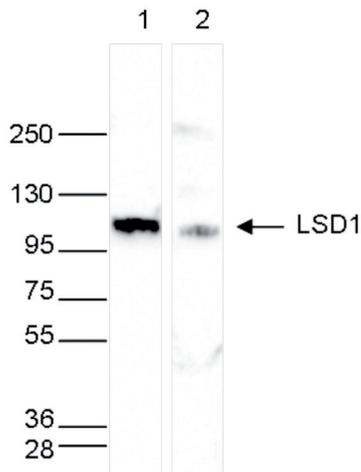


Figure 2. Western blot analysis using the Diagenode antibody directed against LSD1

Whole cell extracts (40 µg) from HeLa cells transfected with LSD1 siRNA (lane 2) and from an untransfected control (lane 1) were analysed by Western blot using the Diagenode antibody against LSD1 (cat. No. C15310028) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.