

PRODUCT NAME		
LSD1 polyclonal antibody		
Other names: BHC110, AOF2, EC1, KDM1		
Cat. No. <b>C15410028</b> (pAb-028-050)	Type: Polyclonal <b>ChIP-grade</b>	Size: 50 µg/ 25 µl
Lot #: 001	Source: Rabbit	Concentration: 2.0 µg/µl

**Description:** Polyclonal antibody raised in rabbit against human LSD1 (Lysine-specific demethylase 1), using the full length recombinant protein.

**Specificity:** Human: positive  
Other species: not tested

Applications	Suggested dilution	References
Western blotting	1:1,000	Fig1
ChIP	3 µg per ChIP	Ref 1

**Purity:** Protein G purified polyclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300.

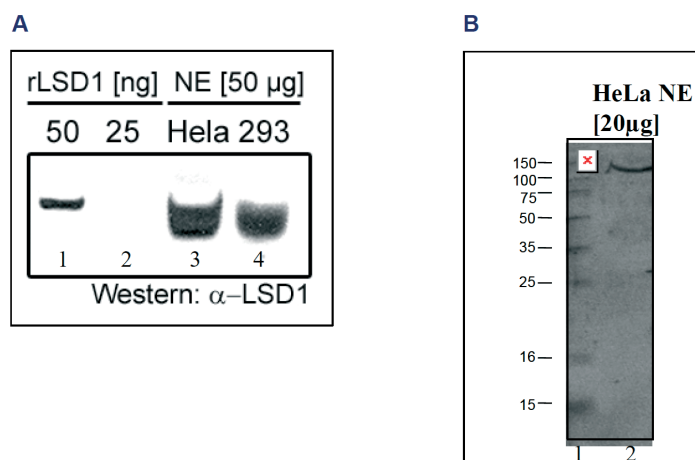
**Storage:** Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

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

**Last data sheet update:** April 9, 2010

#### 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.



**Figure 1**

**Western blot analysis using the Diagenode antibody directed against LSD1**

Western blot was performed on nuclear extracts from HeLa cells (20 µg) using the Diagenode antibody against LSD1 (Cat. No.:pAb-028-050) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The molecular weight marker (in kDa) is shown on the left; the location of the protein of interest is indicated on the right.