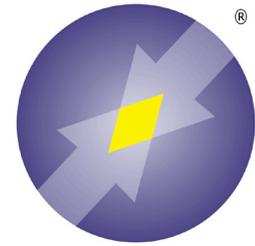


AbX™

## LSD1 Rabbit pAb

Catalog Number A003-100UL



ARBOR  
ASSAYS

### FEATURES

- Reacts with mammalian Lysine Specific Demethylase 1
- Supplied as a PBS solution
- Applications include Western blotting and Immunoprecipitation

### INTRODUCTION

There are two known classes of histone demethylases (HDM), differing in their reaction chemistry, coenzyme use and reaction products. Flavin-dependent HDMs act only on mono- and dimethylated lysines and produce hydrogen peroxide, whereas the *Jumonji*-containing HDMs are iron-dependent enzymes that can act on mono-, di- and trimethylated lysine side chains, and also on methylated Arginine residues. Both classes also produce formaldehyde as a product of the demethylation reaction. LSD1 (Lysine-specific demethylase 1) is a flavin-dependent amine oxidase and was identified as a subunit of different complexes and has been shown to be involved in transcriptional repression of genes through the demethylation of monomethyl Lys<sup>4</sup> and dimethyl Lys<sup>4</sup> Histone H3. LSD1 is also involved in transcriptional activation through the demethylation of monomethyl Lys<sup>9</sup> and dimethyl Lys<sup>9</sup> Histone H3. It is also able to demethylate both mono- and dimethyl Lys<sup>370</sup> in the regulatory domain of the tumor suppressor p53.

### FORM

AbX™ LSD1 Rabbit Antibody is produced as diluted rabbit serum.

### IMMUNOGEN

Full length recombinant Lysine Specific Demethylase 1

### BUFFER COMPOSITION

Phosphate Buffered Saline at pH 7.2 containing 0.1% Tween 20 and 0.09% Kathon preservative

### STORAGE

Short Term: 4°C. Extended: Aliquot and freeze at -20°C

### USES

Western blotting and Immunoprecipitation

### SUGGESTED DILUTION

Western blotting, 1:50-1:200

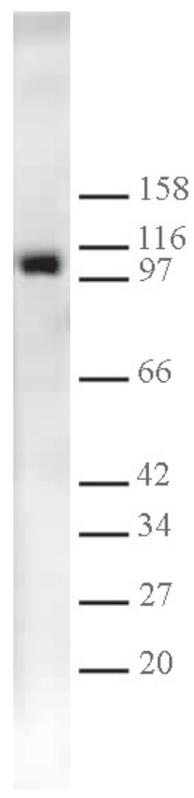
**FOR RESEARCH USE ONLY**

info@ArborAssays.com • www.ArborAssays.com

**WESTERN BLOTTING**

Detection of LSD1 in nuclear extract of 20 µg of HeLa cells probed with A003-100UL at a 1:200 dilution.

MW of LSD1 is 110kDa.



**Related Products**

DetectX™ Histone Demethylase Fluorescent Activity Kit  
Catalog Number K010-F1

AbX™ Antibodies to Unmodified, Monomethylated, Dimethylated and  
Trimethylated Lysine<sup>4</sup> Histone H3  
Catalog Numbers A004, A005, A006, and A007