

## LSD1 / KDM1A antibody (pAb)

**Catalog Nos:** 39186, 39187

**RRID:** AB\_2793175

**Isotype:** Serum

**Application(s):** ChIP, IP, WB

**Reactivity:** Human, Mouse, Wide Range Predicted

**Volume:** 100 µl

**Purification:** None

**Host:** Rabbit

**Molecular Weight:** 110 kDa

**Background:** LSD1 (Lysine-specific demethylase 1, also designated KDM1A) was identified as a subunit of different complexes (such as CTBP, CoREST, NuRD or BRAF35). LSD1 is involved in transcriptional repression of genes through the demethylation of Histone H3K4me1 and Histone H3K4me2, a methylation site associated with transcriptional activation. LSD1 is also involved in transcriptional activation through the demethylation of Histone H3K9me1 and Histone H3K9me2, a modification often associated with transcriptional repression. LSD1 is also able to demethylate Lys370me1 and Lys370me2 in the regulatory domain of the tumor suppressor p53. LSD1 is a flavin-dependent amine oxidase.

**Immunogen:** This antibody was raised against full-length recombinant human LSD1 / KDM1A protein.

**Buffer:** Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

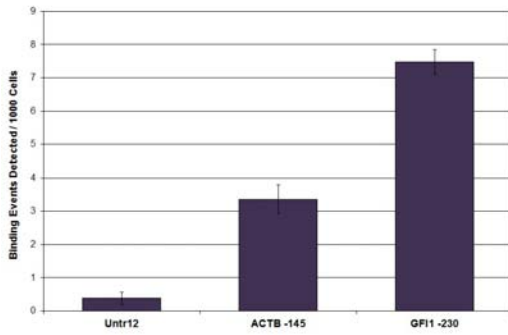
ChIP: 10 µl per ChIP

WB\*: 1:500 - 1:2,000 dilution

\*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

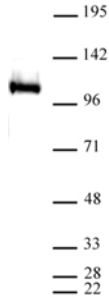
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



**LSD1 / KDM1A antibody (pAb) tested by ChIP.**

Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 30 µg of chromatin from a human small lung cancer cell line (treated with the LSD1 inhibitor, ORY-1001, 1nM for 96 hours) and 5 µl of LSD1 / KDM1A antibody. ChIP DNA was used in qPCR with the negative control primer pairs or gene-specific primer pairs as indicated. Data are presented as Binding Events Detected per 1000 Cells using Active Motifs Epigenetic Services normalization scheme which accounts for primer efficiency and the amount of chromatin used in the ChIP reaction.



**LSD1 / KDM1A antibody (pAb) tested by Western blot**

Nuclear extract of HeLa cells (30 µg per lane) probed with LSD1 / KDM1A antibody (1:2,000 dilution).