

| | | | |
|------------------------------------------|-----------------------------|-------------------------------|-------------------|
| PRODUCT NAME | | | |
| synthetic peptide: inner part of LSD1 | | | |
| (Lysine-specific demethylase 1) | | | |
| Cat. No: sp-067-050 | Format: 50 µg/ 50 µl | Concentration: 1 µg/µl | Lot #: 001 |

Description: Synthetic peptide derived from the inner part of the human LSD1 protein.

This peptide was used as immunogen to raise the Diagenode ChIP-grade antibody directed against LSD1 (cat# pAb-067-050). The same peptide can be used to block this antibody in negative control experiments.

For more information about LSD1: see overview in the Technical Data Sheet of our ChIP-grade antibody directed against LSD1 (cat# pAb-067-050).

Format: In solution in 10 mM HEPES, 2 mM EDTA, 0.05% azide, pH 7.5.

Storage: For long storage, aliquot and store at -20°C. Avoid multiple freeze-thaw cycles.

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

Availability date: November 12, 2007

Last data sheet update: November 12, 2007

Lot #: 001/ day of the synthesis: April 11, 2006/ aliquoting day: November 12, 2007