

<b>PRODUCT NAME</b>			
synthetic peptide: human LSD1			
<b>(Lysine-specific demethylase 1)</b>			
<b>Cat. No:</b> sp-125-050	<b>Format:</b> 50 µg/ 50 µl	<b>Concentration:</b> 1 µg/µl	<b>Lot #:</b> 001

**Description:** Synthetic peptide containing a sequence from the N-terminal part of human LSD1. This peptide was used as immunogen to raise the Diagenode antibody directed against LSD1 (cat# pAb-125-050). 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 the corresponding Diagenode antibody (cat# pAb-125-050).

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

**Storage:** Aliquot upon receipt 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:** July 31, 2009

**Last data sheet update:** July 31, 2009

**Lot #:** 001/ Synthesis date: April 14, 2006/ aliquoting date: July 31, 2009