

Synthetic peptide: human LSD1

(Lysine-specific demethylase 1)

Cat. No. C16000125

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 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. No. C15410125). 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. No. C15410125).

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