

LSD1 (His)

CATALOG NO.: PDM-11-350

LOT NO.:

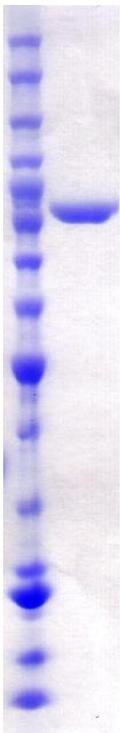
DESCRIPTION: Human recombinant LSD1 (residues 151-852; Genbank Accession # NM_015013.2; MW = 80.3 kDa) expressed with an N-terminal His-tag in *E. coli*.

PURITY: >95% by SDS-PAGE

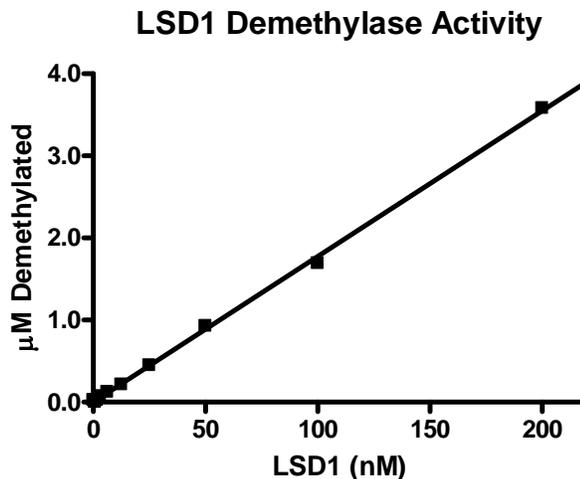
ASSAY CONDITIONS: RBC's LSD1 displays demethyltransferase activity with H3(1-21)K4me2. Reactions were performed with 10 μ M peptide at room temperature in 50mM Tris/HCl, pH 7.5, 0.05% CHAPs, 1% DMSO. H₂O₂ production was monitored by the addition of HRP/Amplex Red mixture to a final concentration of 0.1U/mL and 10 μ M, respectively. The resulting fluorescence signal (Ex. 535 nm/Em. 590 nm) was read every 5 minutes until signal stabilized. (See figure, below.)

SUPPLIED AS: $_ \mu$ g/ μ L in 50 mM Tris/HCl, pH 7.5, 500 mM NaCl, 10% glycerol as determined by OD₂₈₀

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 μ l) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 2 μ g of RBC LSD1 (His). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, & 10 kDa.



LSD1 Demethylase Activity Assay. H3(1-21)K4me2 peptide demethylation was monitored using detection of H₂O₂ production. The 10 μ L reaction contained 10 μ M peptide and a variable concentration of LSD1. H₂O₂ production was detected by the addition of a HRP/Amplex Red detection mixture after 60 minutes. Fluorescence emission (590 nm) was read on an EnVision reader (Perkin Elmer).

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.