

Recombinant LSD1 / KDM1A protein

Catalog No: 31334

Expressed In: Sf9 cells

Quantity: 50 µg

Concentration: 0.5 µg/µl

Source: Human

Buffer Contents: 50 µg recombinant LSD1 / KDM1A protein (greater than 95% purity) expressed in Sf9 cells supplied at a concentration of 0.5 µg/µl in 20 mM Tris-HCl (pH7.3), 300 mM KCl, 0.2 mM EDTA and 20% Glycerol

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 H3 monomethyl Lys4 and Histone H3 dimethyl Lys4, a methylation site associated with transcriptional activation. **LSD1** is also involved in transcriptional activation through the demethylation of Histone H3 monomethyl Lys9 and Histone H3 dimethyl Lys9, 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.

Protein Details: LSD1 (KDM1A, AOF2) is a histone demethylase that is specific to histone H3 dimethylated at lysine 4 (H3K4me2). Methylation of H3K4 is a marker of transcriptional activation. Recombinant full-length human LSD1 (accession number O60341) was expressed with an N-terminal FLAG tag. The molecular weight of the protein is ~95 kDa.

Application Notes: Recombinant LSD1 / KDM1A is suitable for use in histone demethylation assays. A recommended starting point for the demethylation assay is 20 -100 ng of recombinant protein per reaction.

Storage and Guarantee: This product is guaranteed for 6 months from date of receipt.

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