## LSD1 (human), (recombinant)

Highly active

The mono, di- and trimethylation of particular lysine residues in histone tails (e.g. histone H3 lysine-4 (H3K4), H3K9, H3K27, H3K36, H4K20) are implicated, along with a variety of other post-translational modifications (e.g. lysine acetylation) in the transmission of heritable epigenetic information and the control of chromatin structure and DNA transcription. LSD1 catalyzes the oxidative demethylation of mono- and dimethylated H3K4, producing hydrogen peroxide and formaldehyde in the process. H3K4 methylation is considered a transcription activating chromatin mark and *in vivo* LSD1 is frequently found in association with the transcriptional corepressor protein CoREST and HDACs 1 or 2. LSD1 is inhibited by a number of established monoamine oxidase inhibitor drugs. That and the fact that its expression is elevated in a number of cancers may make it a promising target for drug development.

Citations: 4

View Online »

**Ordering Information** 

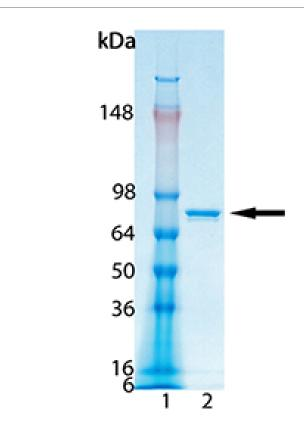
Order Online »

BML-SE544-0050

50µg

Manuals, SDS & CofA

**View Online »** 



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 0.5  $\mu g$  of Purified LSD1 (human) (rec.) Prod. No. BML-SE544.

## **Handling & Storage**

**Use/Stability**The enzyme is stable on ice for the time typically required to set up an experiment (30-

60 min.), but may lose activity with prolonged storage on ice. It is recommended that thawing and dilution of the enzyme be done within as short a time as possible before start of the assay. The remaining, unused, undiluted enzyme should be refrozen quickly by, for example, snap freezing in a dry/ice ethanol bath or liquid nitrogen. Freezing and

storage of diluted enzyme is not recommended.

**Handling** Do not freeze/thaw. After opening, prepare aliquots and store at -80°C.

Long Term Storage -80°C

Shipping Dry Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Activity** Highly active in a peroxidase-coupled assay with the

Histone H3 Dimethyl Lysine-4 Peptide (Prod. No. BML-

P256).

Alternative Name KDM1, Lysine-specific histone demethylase 1, BHC110,

AOF2, KIAA0601

Application Notes Useful tool to study LSD1 kinetics and inhibitor sensitivity

and also the effect of the enzyme in transcriptional regulation, cell cycle progression, and oncogenesis.

Formulation Liquid. In 8.1mM  $Na_2HPO_4$ , 1.5mM  $KH_2PO_4$ , pH 7.2,

138mM sodium chloride, 2.7mM KCl and 40% v/v glycerol.

MW 78 kDa

Purity ≥90% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Sequence N-terminal truncation of LSD1 from human cDNA.

Sequence is identical to Genbank accession NM015013

(aa 151-852).

Source Produced in E. coli.

**Specific Activity** 

≥20 U/µg. One U=1 pmol/min at 30°C, 100 µM Histone H3 Dimethyl Lysine-4 Peptide (Prod. No. BML-P256).

**UniProt ID** 

O60341

Last modified: May 29, 2024



ENZO LIFE SCIENCES, European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com eu@enzolifesciences.com

Phone: 800.942.0430

info-

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com France Phone: +33 472 440 655 fr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 uk@enzolifesciences.com