

# PRODUCT INFORMATION

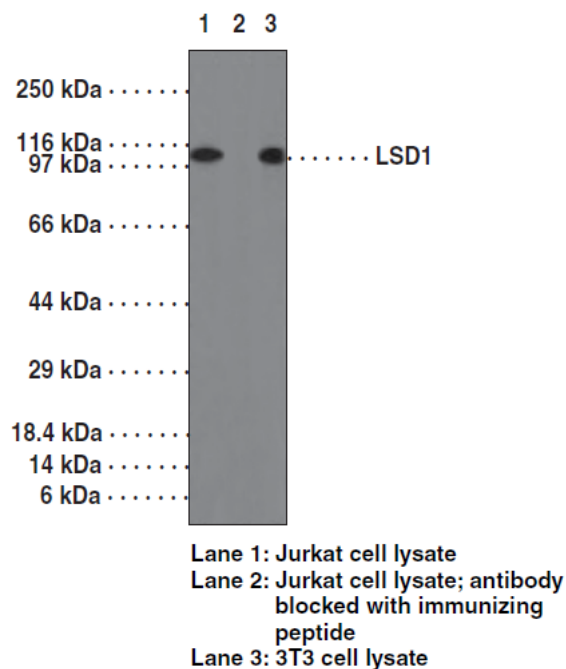
## LSD1 Polyclonal Antibody (aa 450-500)

Item No. 13486

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µg of protein A-purified IgG antibody.
<b>Synonym:</b>	KDM1
<b>Immunogen:</b>	Synthetic peptide from human LSD1 amino acids 450-500
<b>Species Reactivity:</b>	(+) Human, bovine, canine, chimpanzee, equine, mouse, orangutan, porcine; other species not tested
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	200 µl PBS with 0.05% BSA and 0.05% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	Western blot (WB); the recommended starting dilution is 2-3 µg/ml. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/23/2019

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

LSD1, the first known lysine-specific histone demethylase, is an 886 amino acid nuclear protein belonging to flavin monoamine oxidase family. It contains a SWIRM domain, a FAD-binding motif and an amine oxidase domain. This protein is ubiquitously expressed and is a component of several histone deacetylase complexes. LSD1 acts as a component of the CoREST and other transcriptional co-repressor complexes and also plays an important role in silencing neuronal-specific genes in non-neuronal cells. It is also known to demethylate Lys4 of histone H3, a specific tag for epigenetic transcriptional activation. Reports suggest that it plays an important role in stimulating androgen-receptor-dependent transcription converting oxygen to hydrogen peroxide (might use alternative electron acceptors). Along with nuclear FHL2 it serves as a novel biomarker predictive for prostate cancer.

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM