

# Product Information

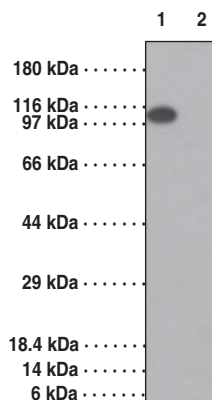


## LSD1 Polyclonal Antibody (aa 100-150)

Item No. 13554

- Synonym:** Amine Oxidase (flavin containing) Domain 2  
**Contents:** This vial contains 100 µg of peptide affinity-purified IgG in 200 µl PBS containing 0.2% gelatin and 0.05% sodium azide.  
**Antigen:** Peptide corresponding to a portion of human LSD1 amino acids 100-150  
**Host:** Rabbit  
**Cross Reactivity:** (+) Canine, human, mouse, rat, Rhesus monkey, and zebrafish LSD1  
**Storage:** ≤6 months at 4°C; ≥6 months at -20°C  
**Applications:** The recommended starting dilution for western blot is 0.5-2 µg/ml.

LSD1 is a flavin-containing nuclear protein having a N-terminal SWIRM domain, a C-terminal FAD-binding motif, and an amine oxidase domain. It functions as a transcriptional corepressor and catalyzes the flavin-dependent demethylation of Lys4 of histone 3 resulting in the formation of methyl-free lysine and release of formaldehyde. LSD1 has been found to be typically associated with CoREST and histone deacetylases 1 and 4 and participates in the silencing of endogenous neuron-specific genes.



Lane 1: Human brain homogenate  
Lane 2: Human brain homogenate-probed with antibody incubated with immunizing peptide

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/13554](http://www.caymanchem.com/catalog/13554)

### Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
[www.caymanchem.com](http://www.caymanchem.com)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 01/29/2013