

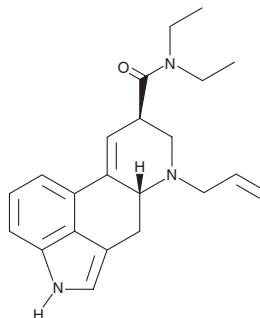
PRODUCT INFORMATION



AL-LAD (solution)

Item No. 30442

CAS Registry No.: 65527-61-9
Formal Name: (8 β)-9,10-didehydro-N,N-diethyl-6-(2-propen-1-yl)-ergoline-8-carboxamide
Synonym: 6-Allyl-nor LSD
MF: C₂₂H₂₇N₃O
FW: 349.5
Supplied as: A 100 μ g/ml solution in acetonitrile
Storage: -20°C
Stability: ≥ 1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

AL-LAD (solution) (Item No. 30442) is an analytical reference material categorized as a lysergamide.¹ AL-LAD substitutes for d-LSD tartrate (Item No. 28289) in a drug discrimination assay in rats.² This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

References

1. Brandt, S.D., Kavanagh, P.V., Westphal, F., *et al.* Return of the lysergamides. Part II: Analytical and behavioural characterization of N⁶-allyl-6-norlysergic acid diethylamide (AL-LAD) and (2'S,4'S)-lysergic acid 2,4-dimethylazetidide (LSZ). *Drug Test Anal.* **9**(1), 38-50 (2017).
2. Hoffman, A.J. and Nichols, D.E. Synthesis and LSD-like discriminative stimulus properties in a series of N(6)-alkyl norlysergic acid N,N-diethylamide derivatives. *J. Med. Chem.* **28**(9), 1252-1255 (1985).

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, 10/05/2020

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