

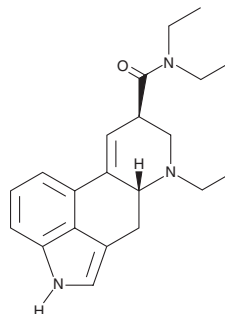
PRODUCT INFORMATION



ETH-LAD (solution)

Item No. 41271

CAS Registry No.: 65527-62-0
Formal Name: 9,10-didehydro-N,N,6-triethyl-ergoline-8 β -carboxamide
Synonyms: N⁶-ethyl-nor-LAD, N⁶-ethyl-nor-Lysergic Acid diethylamide
MF: C₂₁H₂₇N₃O
FW: 337.5
Purity: \geq 95%
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: \geq 2 years



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

Description

ETH-LAD (solution) (Item No. 41271) is an analytical reference standard categorized as a lysergamide.¹ ETH-LAD substitutes for LSD (D-tartrate) (Item No. 28289) in a two-lever drug discrimination test in rats.² This product is intended for research and forensic applications.

References

1. Brandt, S.D., Kavanagh, P.V., Westphal, F., *et al.* Return of the lysergamides. Part III: Analytical characterization of N⁶-ethyl-6-norlysergic acid diethylamide (ETH-LAD) and 1-propionyl ETH-LAD (1P-ETH-LAD). *Drug Test Anal.* **9(10)**, 1641-1649 (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, 02/28/2025

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