

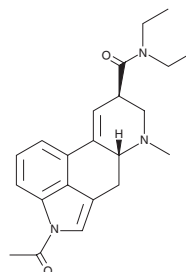
PRODUCT INFORMATION



ALD-52 (solution)

Item No. 27818

CAS Registry No.: 3270-02-8
Formal Name: 1-acetyl-9,10-didehydro-N,N-diethyl-6-methyl-ergoline-8 β -carboxamide
Synonyms: 1-Acetyllysergic Acid diethylamide, 1-acetyl LSD
MF: C₂₂H₂₇N₃O₂
FW: 365.5
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

ALD-52 (solution) (Item No. 27818) is an analytical reference material categorized as a lysergamide.¹ ALD-52 has psychotomimetic properties that are equipotent to those of LSD.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 I: Analytical and behavioural characterization of 1-propionyl-D-lysergic acid diethylamide (1P-LSD). *Drug Test Anal.* **8**(9), 891-902 (2016).
2. Isbell, H., Miner, E.J., and Logan, C.R. Relationships of psychotomimetic to anti-serotonin potencies of congeners of lysergic acid diethylamide (LSD-25). *Psychopharmacologia* **1**, 20-28 (1959).

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, 07/10/2019

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