

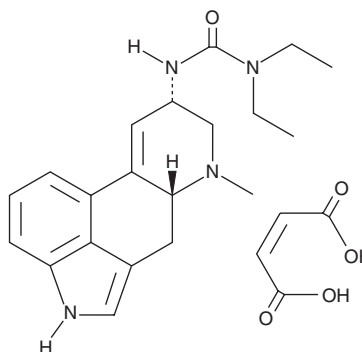
PRODUCT INFORMATION



Lisuride (maleate)

Item No. 36140

CAS Registry No.: 19875-60-6
Formal Name: N'-[(8 α)-9,10-didehydro-6-methylergolin-8-yl]-N,N-diethyl-urea, 2Z-butenedioate
Synonym: Lysenyl
MF: C₂₀H₂₆N₄O • C₄H₄O₄
FW: 454.5
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Light sensitive



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

Description

Lisuride (maleate) (Item No. 36140) is an analytical reference standard categorized as a lysergamide. Lisuride does not induce the head-twitch response (HTR) in mice, indicating a lack of hallucinogenic potential.¹ This product is intended for research and forensic applications.

Reference

1. Halberstadt, A.L. and Geyer, M.A. Characterization of the head-twitch response induced by hallucinogens in mice: Detection of the behavior based on the dynamics of head movement. *Psychopharmacology (Berl)* 227(4), 727-739 (2013).

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, 01/14/2026

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