

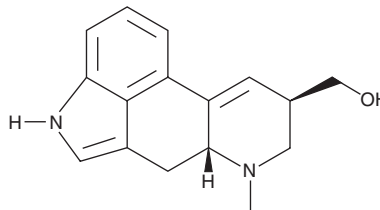
PRODUCT INFORMATION



Lysergol

Item No. 11717

CAS Registry No.: 602-85-7
Formal Name: 9,10-didehydro-6-methyl-ergoline-8 β -methanol
Synonyms: (+)-Lysergol, NSC 196867
MF: C₁₆H₁₈N₂O
FW: 254.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 227, 242, 311 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Fungal



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

Laboratory Procedures

Lysergol is supplied as a crystalline solid. A stock solution may be made by dissolving the lysergol in the solvent of choice, which should be purged with an inert gas. Lysergol is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of lysergol in these solvents is approximately 20 mg/ml.

Description

Lysergol is an ergot alkaloid that has been found in *S. sorghi* and an agonist of the serotonin (5-HT) receptor subtypes 5-HT_{1A}, 5-HT_{2A}, and 5-HT_{2C} (EC₅₀s = 73, 1.6, and 6.6 nM, respectively, in a FLIPR assay).^{1,2} It induces contractions in isolated rat tail arteries (EC₅₀ = 436.5 nM).² Lysergol (10 μ g/ml) also enhances the antibacterial activity of the quinolone antibiotic nalidixic acid (Item No. 19807) against susceptible and drug-resistant *E. coli*.³

References

1. Yuan, H., Guo, Z., and Luo, T. Synthesis of (+)-lysergol and its analogues to assess serotonin receptor activity. *Org. Lett.* **19(3)**, 624-627 (2017).
2. Pertz, H. Naturally occurring clavines: Antagonism/partial agonism at 5-HT_{2A} receptors and antagonism at α_1 -adrenoceptors in blood vessels. *Planta Med.* **62(5)**, 387-392 (1996).
3. Maurya, A., Dwivedi, G.R., Darokar, M.P., et al. Antibacterial and synergy of clavine alkaloid lysergol and its derivatives against nalidixic acid-resistant *Escherichia coli*. *Chem. Biol. Drug Des.* **81(4)**, 484-490 (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, 07/24/2023

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