

PRODUCT INFORMATION



(-)-Himgaline

Item No. 29990

CAS Registry No.: 16585-08-3
Formal Name: (1R,4S,5S,5aS,7R,7aR,11aS,11bS,12S,13aR,15R)-hexadecahydro-4-methyl-12,1,5a-ethanylylidene-5aH-naphtho[1,2-c]quinolizine-7,15-diol

MF: C₂₀H₃₁NO₂

FW: 317.5

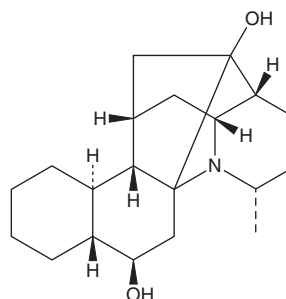
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Galbulimima* sp.



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

Laboratory Procedures

(-)-Himgaline is supplied as a solid. A stock solution may be made by dissolving the (-)-himgaline in the solvent of choice, which should be purged with an inert gas. (-)-Himgaline is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

(-)-Himgaline is an alkaloid that has been found in *G. belgraveana*.¹

Reference

1. Shah, U., Chackalamannil, S., Ganguly, A.K., *et al.* Total synthesis of (-)-himgaline. *J. Am. Chem. Soc.* **128**(39), 12654-12655 (2006).

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, 10/25/2022

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