

PRODUCT INFORMATION



Himandridine

Item No. 30055

CAS Registry No.: 15521-76-3
Formal Name: (1S,2R,4aS,4bS,5R,6S,7aR,9S,11R,11aR,12R,15R)-11-(benzoyloxy)-1,2,3,4,4a,6,7,7a,8,9,11,11a-dodecahydro-2,9-dihydroxy-1-methoxy-15-methyl-5,6,9-[1]Butanyl[4]ylidene-5H-benz[6,7]indeno[7a,1-b]pyrrole-10-carboxylic acid, methyl ester

Synonym: NSC 23972

MF: C₃₀H₃₇NO₇

FW: 523.6

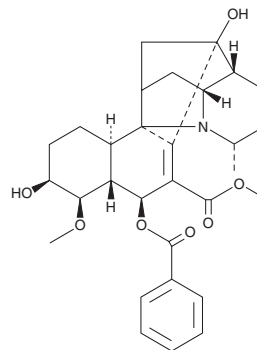
Purity: ≥95%

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

Himandridine is supplied as a solid. A stock solution may be made by dissolving the himandridine in the solvent of choice, which should be purged with an inert gas. Himandridine is soluble in ethanol, methanol, DMSO, and dimethylformamide.

Description

Himandridine is an alkaloid that has been found in *Galbulimima*.¹

Reference

1. Mander, L.N., Ritchie, E., and Taylor, W.C. The chemical constituents of *Galbulimima* species. VI. The structure of himandridine. *Aust. J. Chem.* **20(5)**, 981-1019 (1967).

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, 10/27/2022

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