

PRODUCT INFORMATION



Emindole SB

Item No. 44964

CAS Registry No.: 112900-04-6
Formal Name: 1,2,3S,4S,4aR,5,6,6aS,7,12,12bS,12cS-dodecahydro-4,12b,12c-trimethyl-4-(4-methyl-3-penten-1-yl)-benz[6,7]indeno[1,2-b]indol-3-ol

MF: C₂₈H₃₉NO

FW: 405.6

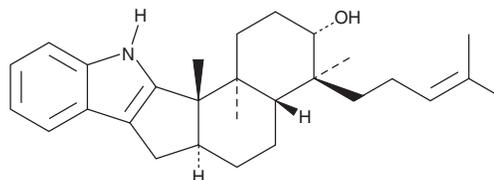
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungi/*Aspergillus* sp.



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

Laboratory Procedures

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

Description

Emindole SB is an indoloditerpene that has been found in *D. cejpaii* and has diverse biological activities.¹ It is a cannabinoid 1 (CB₁) and CB₂ receptor antagonist (K_is = 7.05 and 2.24 μM, respectively). Emindole SB inhibits protein tyrosine phosphatase 1B (PTP1B) and T cell protein tyrosine phosphatase (TCPTP; IC₅₀s = 1.6 and 2.4 μM, respectively).²

References

1. Harms, H., Rempel, V., Kehraus, S., *et al.* Indoloditerpenes from a marine-derived fungal strain of *Dichotomomyces cejpaii* with antagonistic activity at GPR18 and cannabinoid receptors. *J. Nat. Prod.* **77(3)**, 673-677 (2014).
2. Kong, F.-D., Fan, P., Zhou, L.-M., *et al.* Peneperenes A-D, four indole terpenoids with potent protein tyrosine phosphatase inhibitory activity from the marine-derived fungus *Penicillium* sp. KFD28. *Org. Lett.* **21(12)**, 4864-4867 (2019).

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, 01/23/2026

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