

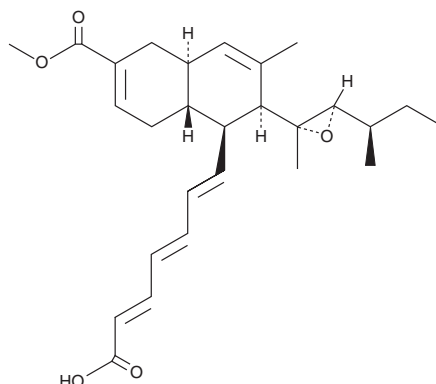
PRODUCT INFORMATION



Cladobotric Acid H

Item No. 45500

CAS Registry No.: 2900278-39-7
Formal Name: (4aR,5S,6S,8aR)-5-[(1E,3E,5E)-6-carboxy-1,3,5-hexatrien-1-yl]-1,4,4a,5,6,8a-hexahydro-7-methyl-6-[(2R,3R)-2-methyl-3-[(1R)-1-methylpropyl]-2-oxiranyl]-2-naphthalenecarboxylic acid, 2-methyl ester
MF: C₂₇H₃₆O₅
FW: 440.6
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Cladobotryum* sp.



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

Laboratory Procedures

Cladobotric acid H is supplied as a crystalline solid. A stock solution may be made by dissolving the cladobotric acid H in the solvent of choice, which should be purged with an inert gas. Cladobotric acid H is slightly soluble (0.1-1 mg/ml) in methanol.

Description

Cladobotric acid H is a fungal metabolite that has been found in *H. pseudocorticiicola* with antibacterial and antifungal activities.^{1,2} It is active against methicillin-susceptible, methicillin-resistant and vancomycin-intermediate, and heterogeneous vancomycin-intermediate strains of *S. aureus* (MICs = 64 µg/ml for all).¹ Cladobotric acid H is also active against both azole-sensitive and azole-resistant strains of *A. fumigatus* (MICs = 32 µg/ml) and *C. auris* (MICs = 16-32 µg/ml).²

References

1. Dao, T.T., Williams, K., de Mattos-Shiple, K.M.J., et al. Cladobotric acids: Metabolites from cultures of *Cladobotryum* sp., semisynthetic analogues and antibacterial activity. *J. Nat. Prod.* **85**(3), 572-580 (2022).
2. Watanabe, Y., Takahashi, S., Ito, S., et al. Hakuhybotrol, a polyketide produced by *Hypomyces pseudocorticiicola*, characterized with the assistance of 3D ED/MicroED. *Org. Biomol. Chem.* **21**(11), 2320-2330 (2023).

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, 04/02/2026

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