

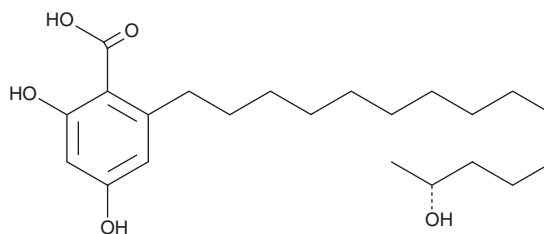
PRODUCT INFORMATION



Phanerosporic Acid

Item No. 23879

CAS Registry No.: 124709-28-0
Formal Name: 2,4-dihydroxy-6-[(14R)-14-hydroxypentadecyl]-benzoic acid
Synonym: (-)-Phanerosporic Acid
MF: C₂₂H₃₆O₅
FW: 380.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Phanerochaete* sp.



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

Laboratory Procedures

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

Description

Phanerosporic acid is a fungal secondary metabolite originally isolated from *P. chrysosporum* that has antibacterial activity against *B. cereus*, *B. subtilis*, and *E. coli* and antifungal activity against *S. cerevisiae*, *A. niger*, *O. ulmi*, *U. maydis*, *C. cucumerinum*, *C. cladosporioides*, and *B. cinerea*.¹ It has been used in the synthesis of macrolide derivatives.

Reference

1. Arnone, A., Assante, G., Nasini, G., *et al.* Phanerosporic acid, a β -resorcyolate obtained from *Phanerochaete chrysosporium*. *Phytochemistry* **28(10)**, 2803-2806 (1989).

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, 08/23/2023

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