

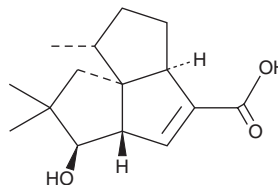
PRODUCT INFORMATION



Pentalenic Acid

Item No. 33532

CAS Registry No.: 69394-19-0
Formal Name: (1R,3aR,5aS,6R,8aS)-1,2,3,3a,5a,6,7,8-octahydro-6-hydroxy-1,7,7-trimethyl-cyclopenta[c]pentalene-4-carboxylic acid
MF: $C_{15}H_{22}O_3$
FW: 250.3
Purity: $\geq 95\%$
Supplied as: An oil
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Bacterium/*Streptomyces UC5319*



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

Laboratory Procedures

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

Description

Pentalenic acid is a sesquiterpenoid that has been found in *S. avermitilis*.¹ It is a precursor in the biosynthesis of the antibiotic pentalenolactone.

Reference

1. Takamatsu, S., Xu, L.-H., Fushinobu, S., *et al.* Pentalenic acid is a shunt metabolite in the biosynthesis of the pentalenolactone family of metabolites: Hydroxylation of 1-deoxypentalenic acid mediated by CYP105D7 (SAV_7469) of *Streptomyces avermitilis*. *J. Antibiot.* **64**(1), 65-71 (2011).

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, 02/19/2025

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