

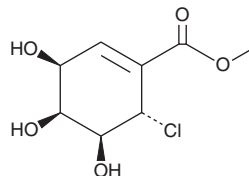
PRODUCT INFORMATION



Pericosine A

Item No. 29633

CAS Registry No.: 200335-68-8
Formal Name: (3S,4S,5S,6S)-6-chloro-3,4,5-trihydroxy-1-cyclohexene-1-carboxylic acid, methyl ester
Synonym: (+)-Pericosine A
MF: C₈H₁₁ClO₅
FW: 222.6
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

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

Description

Pericosine A is a fungal metabolite that has been found in *P. byssoides* and has anticancer activity.¹ It inhibits the growth of a variety of cancer cells, including breast, colon, lung, ovary, stomach, and prostate cell lines (GI₅₀s = 0.05-24.55 μM) and increases survival in a P388 mouse xenograft model when administered at a dose of 25 mg/kg. Pericosine A inhibits EGFR by 40 to 70% when used at a concentration of 100 μg/ml. It also reacts with organosulfur compounds in skunk spray to form stable thioethers as odorless products.²

References

1. Yamada, T., Iritani, M., Ohishi, H., *et al.* Pericosines, antitumour metabolites from the sea hare-derived fungus *Periconia byssoides*. Structures and biological activities. *Org. Biomol. Chem.* **5(24)**, 3979-3986 (2007).
2. Du, L., Munteanu, C., King, J.B., *et al.* An electrophilic natural product provides a safe and robust odor neutralization approach to counteract malodorous organosulfur metabolites encountered in skunk spray. *J. Nat. Prod.* **82(7)**, 1989-1999 (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, 08/15/2023

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