

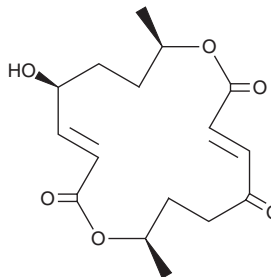
PRODUCT INFORMATION



Dihydropyrenophorin

Item No. 33536

CAS Registry No.: 1073287-13-4
Formal Name: 13S-hydroxy-8R,16R-dimethyl-1,9-dioxacyclohexadeca-3E,11E-diene-2,5,10-trione
Synonym: (-)-Dihydropyrenophorin
MF: C₁₆H₂₂O₆
FW: 310.3
Purity: ≥95%
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

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

Description

Dihydropyrenophorin is a macrolide fungal metabolite and derivative of pyrenophorin (Item No. 33502) originally isolated from *D. avenae* that has herbicidal and phytotoxic activities.¹ It is herbicidal against *S. halepense* when used at a concentration of 0.065 μM. Dihydropyrenophorin is phytotoxic to a variety of agricultural crops, including barley, sorghum, wheat, and oats, as well as grasses, including goosegrass and Bermuda grass, in leaf puncture assays.

Reference

1. Sugawara, F. and Strobel, G.A. (-)-Dihydropyrenophorin, a novel and selective phytotoxin produced by *Drechslera avenae*. *Plant Sci.* **43(1)**, 1-5 (1986).

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, 08/25/2023

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