

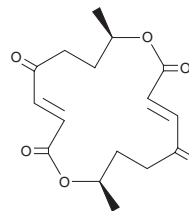
PRODUCT INFORMATION



Pyrenophorin

Item No. 33502

CAS Registry No.: 5739-85-5
Formal Name: (3E,8R,11E,16R)-8,16-dimethyl-1,9-dioxacyclohexadeca-3,11-diene-2,5,10,13-tetrone
Synonym: (-)-Pyrenophorin
MF: C₁₆H₂₀O₆
FW: 308.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Description

Pyrenophorin is a fungal metabolite that has been found in *D. avenae* and has phytotoxic activity.¹ It inhibits seed germination and reduces root growth of the graminaceous plants *A. sterilis*, *A. fatua*, *A. sativa*, *T. aestivum*, and *H. vulgare* when used at a concentration of 60 μM. Pyrenophorin (320 μM) reduces chlorophyll levels in leaf sections from seedlings of the same plants.

Reference

1. Kastanias, M.A. and Chrysayi-Tokousbalides, M. Comparative phytotoxicity of pyrenophorin and pyrenophorol isolated from *Drechslera avenae* pathotype. 1999 Brighton crop protection conference: Weeds 2, 581-582 (1999).

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, 11/15/2022

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