

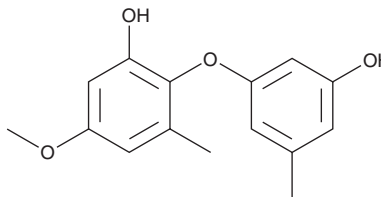
PRODUCT INFORMATION



Cyperine

Item No. 29937

CAS Registry No.: 33716-82-4
Formal Name: 2-(3-hydroxy-5-methylphenoxy)-5-methoxy-3-methyl-phenol
Synonym: Cyperin
MF: C₁₅H₁₆O₄
FW: 260.3
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Zopfia* sp.



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

Laboratory Procedures

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

Description

Cyperine is a phytotoxic fungal metabolite that has been found in *A. cypericola*.¹ It inhibits *A. thaliana* enoyl-acyl carrier protein reductase (ENR; IC₅₀ = 89 μM).² Cyperine induces necrosis in a panel of nine plant species in a leaf bioassay and inhibits root growth in *A. thaliana* seedlings (IC₅₀ = 38.4 μM).^{2,3}

References

1. Stierle, A., Upadhyay, R.S., and Strobel, G. Cyperine, a phytotoxin produced by *Ascochyta cypericola*, a fungal pathogen of *Cyperus rotundus*. *Phytochemistry* **30(7)**, 2191-2192 (1991).
2. Dayan, F.E., Ferreira, D., Wang, Y.-H., et al. A pathogenic fungi diphenyl ether phytotoxin targets plant enoyl (acyl carrier protein) reductase. *Plant Physiol.* **174(3)**, 1062-1071 (2008).
3. Venkatasubbaiah, P., Van Dyke, C.G., and Chilton, W.S. Phytotoxic metabolites of *Phoma sorghina*, a new foliar pathogen of pokeweed. *Mycologia* **84(5)**, 715-723 (1992).

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, 08/16/2023

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