

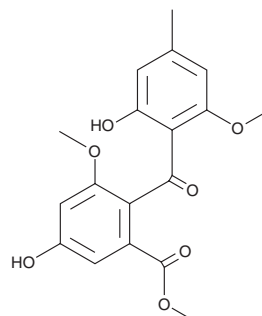
PRODUCT INFORMATION



Monomethylsulochrin

Item No. 29942

CAS Registry No.: 10056-14-1
Formal Name: 5-hydroxy-2-(2-hydroxy-6-methoxy-4-methylbenzoyl)-3-methoxy-benzoic acid, methyl ester
Synonyms: 1,2-Secotrypacidin, Methylsulochrin
MF: C₁₈H₁₈O₇
FW: 346.3
Purity: ≥98%
UV/Vis.: λ_{max}: 284 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungi/*Aspergillus fumigatus*



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

Laboratory Procedures

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

Description

Monomethylsulochrin is a fungal metabolite that has been found in *Rhizoctonia*.¹ It is active against *H. pylori* (MIC = 10 µg/ml).

Reference

1. Ma, Y.M., Li, Y., Liu, J.Y., et al. Anti-*Helicobacter pylori* metabolites from *Rhizoctonia* sp. Cy064, an endophytic fungus in *Cynodon dactylon*. *Fitoterapia* **75**(5), 451-456 (2004).

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, 12/12/2022

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