

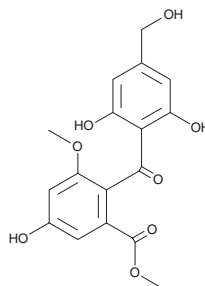
PRODUCT INFORMATION



Hydroxysulochrin

Item No. 29325

CAS Registry No.: 581068-64-6
Formal Name: 2-[2,6-dihydroxy-4-(hydroxymethyl)benzoyl]-5-hydroxy-3-methoxybenzoic acid, methyl ester
MF: $C_{17}H_{16}O_8$
FW: 348.3
Purity: $\geq 70\%$
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

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

Description

Hydroxysulochrin is a fungal metabolite and plant growth regulator that has been found in *Aureobasidium*.¹ It inhibits tea (*C. sinensis*) pollen tube growth by 41%, but does not inhibit lettuce (*L. sativa*) seedling root growth, when used at a concentration of 100 mg/L.

Reference

1. Shimada, A., Shiokawa, C., Kusano, M., *et al.* Hydroxysulochrin, a tea pollen growth inhibitor from the fungus *Aureobasidium* sp. *Biosci. Biotechnol. Biochem.* **67**(2), 442-444 (2003).

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