

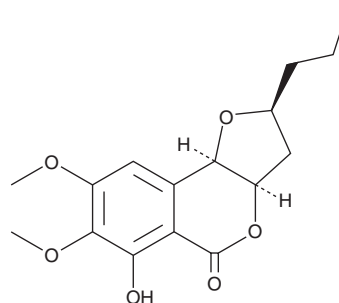
PRODUCT INFORMATION



Monocerin

Item No. 29572

CAS Registry No.: 30270-60-1
Formal Name: (2S,3aR,9bR)-2,3,3a,9b-tetrahydro-6-hydroxy-7,8-dimethoxy-2-propyl-5H-furo[3,2-c][2]benzopyran-5-one
Synonym: (+)-Monocerin
MF: C₁₆H₂₀O₆
FW: 308.3
Purity: ≥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

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

Description

Monocerin is a fungal metabolite that has been found in *F. larvarum* and has diverse biological activities.¹⁻³ It is active against the bacteria *E. coli* and *B. megaterium*, the phytopathogenic fungus *M. violaceum*, and the alga *C. fusca* in an agar diffusion assay when used at a concentration of 50 µg/disc.¹ Monocerin (17.5 µg/ml) induces mortality in adult *C. erythrocephala*.² It reduces root elongation in pre-germinated *S. halepense* seeds when used at a concentration of 33 ppm.³

References

1. Zhang, W., Krohn, K., Draeger, S., *et al.* Bioactive isocoumarins isolated from the endophytic fungus *Microdochium bolleyi*. *J. Nat. Prod.* **71(6)**, 1078-1081 (2008).
2. Claydon, N., Grove, J.F., and Pople, M. Insecticidal secondary metabolic products from the entomogenous fungus *Fusarium larvarum*. *J. Invertebr. Pathol.* **33(3)**, 364-367 (1979).
3. Robeson, D.J. and Strobel, G.A. Monocerin, a phytotoxin from *Exserohilum turcicum* (*Drechslera turcica*). *Agr. Biol. Chem.* **46(11)**, 2681-2683 (1982).

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