

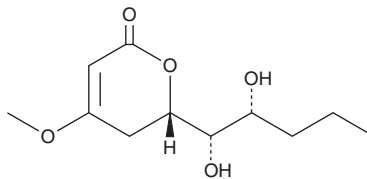
PRODUCT INFORMATION



7-hydroxy Pestalotin

Item No. 26523

CAS Registry No.: 41164-59-4
Formal Name: (6S)-6-[(1S,2R)-1,2-dihydroxypentyl]-5,6-dihydro-4-methoxy-2H-pyran-2-one
MF: $C_{11}H_{18}O_5$
FW: 230.3
Purity: $\geq 99\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*Penicillium decumbens*



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

Laboratory Procedures

7-hydroxy Pestalotin is supplied as a solid. A stock solution may be made by dissolving the 7-hydroxy pestalotin in the solvent of choice, which should be purged with an inert gas. 7-hydroxy Pestalotin is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

7-hydroxy Pestalotin has limited solubility in water. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

7-hydroxy Pestalotin is a fungal metabolite originally isolated from *Penicillium*.¹ It is phytotoxic against evening primrose, prickly sida, johnsongrass, morning glory, lambsquarter, and *A. abla*.²

References

1. McGahren, W.J., Ellestad, G.A., Morton, G.O., *et al.* A new fungal lactone, LL-P880 β , and a new pyrone, LL-P880 γ , from a *Penicillium* sp. *J. Org. Chem.* **38**(20), 3542-3544 (1973).
2. Venkatasubbaiah, P., Van Dyke, C.G., and Chilton, W.S. Phytotoxins, produced by *Pestalotiopsis Oenotherae*, a pathogen of evening primrose. *Phytochemistry* **30**(5), 1471-1474 (1991).

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, 07/07/2025

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