

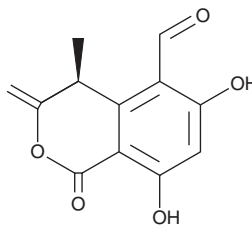
PRODUCT INFORMATION



Ascochin

Item No. 31450

CAS Registry No.: 935699-58-4
Formal Name: (4S)-3,4-dihydro-6,8-dihydroxy-4-methyl-3-methylene-1-oxo-1H-2-benzopyran-5-carboxaldehyde
Synonyms: (4S)-(+)-Ascochin, (S)-(+)-Ascochin
MF: C₁₂H₁₀O₅
FW: 234.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Ascochyta* sp.



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

Laboratory Procedures

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

Description

Ascochin is an isocoumarin derivative and fungal metabolite that has been found in *Ascochyta*.¹ It is active against the bacterium *B. megaterium*, the plant pathogenic fungus *M. violaceum*, and the alga *C. fusca* in an agar diffusion assay when used at a concentration of 0.5 mg/disc.

Reference

1. Kohn, K., Kock, I., Elsässer, B., *et al.* Bioactive natural products from the endophytic fungus *ascochyta* sp. from *melilotus dentatus* – configurational assignment by solid-state CD and TDDFT calculations. *Eur. J. Org. Chem.* 1123-1129 (2007).

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, 11/10/2022

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