

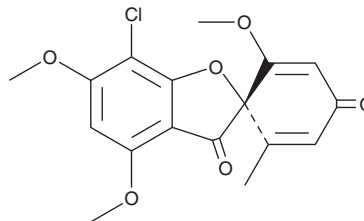
PRODUCT INFORMATION



(-)-dehydro Griseofulvin

Item No. 33117

CAS Registry No.: 3573-90-8
Formal Name: (1'S)-7-chloro-2',4,6-trimethoxy-6'-methyl-spiro[benzofuran-2(3H),1'-[2,5]cyclohexadiene]-3,4'-dione
MF: C₁₇H₁₅ClO₆
FW: 350.8
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

(-)-dehydro Griseofulvin is supplied as a solid. A stock solution may be made by dissolving the (-)-dehydro griseofulvin in the solvent of choice, which should be purged with an inert gas. (-)-dehydro Griseofulvin is slightly soluble in acetone and DMSO.

Description

(-)-dehydro Griseofulvin is a polyketide and derivative of griseofulvin (Item No. 19461) that has been found in *Penicillium*.^{1,2}

References

1. Zhang, H.-M., Ju, C.-X., Li, G., *et al.* Dimeric 1,4-benzoquinone derivatives with cytotoxic activities from the marine-derived fungus *Penicillium* sp. L129. *Mar. Drugs* **17**(7), 383 (2019).
2. Cacho, R.A., Chooi, Y.-H., Zhou, H., *et al.* Complexity generation in fungal polyketide biosynthesis: A spirocycle-forming P450 in the concise pathway to the antifungal drug griseofulvin. *ACS Chem. Biol.* **8**(10), 2322-2330 (2013).

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, 11/14/2022

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