

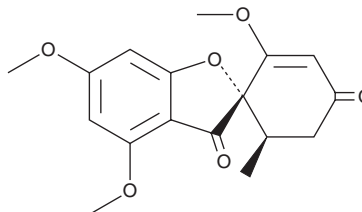
PRODUCT INFORMATION



7-dechloro Griseofulvin

Item No. 33116

CAS Registry No.: 3680-32-8
Formal Name: (1'S,6'R)-2',4,6-trimethoxy-6'-methyl-spiro[benzofuran-2(3H), 1'-[2]cyclohexene]-3,4'-dione
MF: C₁₇H₁₈O₆
FW: 318.3
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

7-dechloro Griseofulvin is supplied as a solid. A stock solution may be made by dissolving the 7-dechloro griseofulvin in the solvent of choice, which should be purged with an inert gas. 7-dechloro Griseofulvin is slightly soluble in acetone, DMSO, and ethyl acetate (warmed).

Description

7-dechloro Griseofulvin is a polyketide and derivative of griseofulvin (Item No. 19461) that has been found in *Penicillium* and has fungicidal and plant growth inhibitory properties.¹⁻³ It is active against the plant pathogenic fungi *B. cinerea*, *C. orbiculare*, *R. solani*, and *S. sclerotiorum* with EC₅₀ values of 40, 81.7, 43.9, and 28.5 µg/ml, respectively.² 7-dechloro Griseofulvin (1-300 mg/L) inhibits hypocotyl and root growth of lettuce seedlings.³

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. Zhao, J.H., Zhang, Y.L., Wang, L.W., *et al.* Bioactive secondary metabolites from *Nigrospora* sp. LLGLM003, an endophytic fungus of the medicinal plant *Moringa oleifera* Lam. *World J. Microbiol. Biotechnol.* **28(5)**, 2107-2112 (2012).
3. Kimura, Y., Shiojima, K., Nakajima, H., *et al.* Structure and biological activity of plant growth regulators produced by *Penicillium* sp. No. 31f. *Biosci. Biotechnol. Biochem.* **56(7)**, 1138-1139 (1992).

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, 12/09/2022

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