

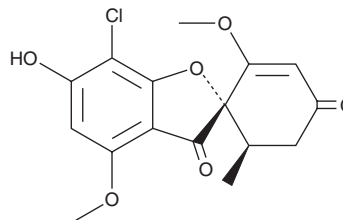
PRODUCT INFORMATION



6-O-Demethyl Griseofulvin

Item No. 33118

CAS Registry No.: 20168-88-1
Formal Name: (1'S,6'R)-7-chloro-6-hydroxy-2',4-dimethoxy-6'-methyl-spiro[benzofuran-2(3H),1'-[2]cyclohexene]-3,4'-dione
Synonym: 6-Demethylgriseofulvin
MF: C₁₆H₁₅ClO₆
FW: 338.7
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Penicillium* sp.



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

Laboratory Procedures

6-O-demethyl Griseofulvin is supplied as a solid. A stock solution may be made by dissolving the 6-O-demethyl griseofulvin in the solvent of choice, which should be purged with an inert gas. 6-O-Demethyl Griseofulvin is soluble in methanol and DMSO.

Description

6-O-demethyl Griseofulvin is a metabolite of the antifungal griseofulvin (Item No. 19461).^{1,2}

References

1. Zia, H., Proveaux, W.J., O'Donnell, J.P., *et al.* Chromatographic analysis of griseofulvin and metabolites in biological fluids. *J. Chromatogr.* **181(1)**, 77-84 (1980).
2. Schäfer-Korting, M., Korting, H.C., and Mutschler, E. Human plasma and skin blister fluid levels of griseofulvin after its repeated administration. *Eur. J. Clin. Pharmacol.* **29(3)**, 351-354 (1985).

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, 08/23/2023

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