

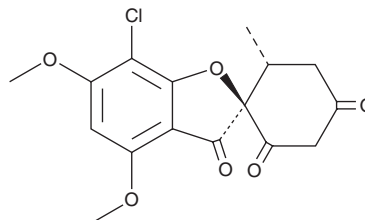
PRODUCT INFORMATION



Griseofulvic Acid

Item No. 33086

CAS Registry No.: 469-54-5
Formal Name: 7-chloro-4,6-dimethoxy-6'R-methyl-spiro[benzofuran-2(3H),1'S-cyclohexane]-2',3,4'-trione
Synonym: (±)-Griseofulvic Acid
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

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

Description

Griseofulvic acid is a metabolite of the antifungal griseofulvin (Item No. 19461).¹ It induces protein aggregation and tubulin polymerization in cell-free assays.²

References

1. Zia, H., O'Donnell, J.P., and Ma, J.K. Identification of griseofulvic acid as a urine metabolite of griseofulvin in humans. *J. Pharm. Sci.* **68**(10), 1335-1336 (1979).
2. Mir, L., Oustrin, M.-L., Lecointe, P., *et al.* Correlation between the in vivo effects of some griseofulvin derivatives and their in vitro interactions with mammalian microtubules. *FEBS Lett.* **88**(2), 259-263 (1978).

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, 09/26/2023

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