

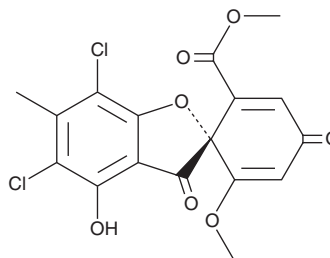
PRODUCT INFORMATION



(+)-Geodin

Item No. 26421

CAS Registry No.: 427-63-4
Formal Name: (1'R)-5,7-dichloro-4-hydroxy-6'-methoxy-6-methyl-3,4'-dioxo-spiro[benzofuran-2(3H),1'-[2,5]cyclohexadiene]-2'-carboxylic acid, methyl ester
MF: C₁₇H₁₂Cl₂O₇
FW: 399.2
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

(+)-Geodin is supplied as a solid. A stock solution may be made by dissolving the (+)-geodin in the solvent of choice, which should be purged with an inert gas. (+)-Geodin is soluble in organic solvents such as methanol, DMSO, chloroform, and ethyl acetate.

Description

(+)-Geodin is a fungal metabolite.^{1,2} It increases the fibrinolytic activity of bovine aortic endothelial cells (BAECs) when used at concentrations ranging from 50 to 150 μM.¹ (+)-Geodin (1-100 μg/ml) also increases 2-deoxyglucose uptake by rat adipocytes.²

References

1. Shinohara, C., Chikanishi, T., Nakashima, S., *et al.* Enhancement of fibrinolytic activity of vascular endothelial cells by chaetoglobosin A, crinipellin B, geodin and triticone B. *J. Antibiot. (Tokyo)* **53**(3), 262-268 (2000).
2. Sato, S., Okusa, N., Ogawa, A., *et al.* Identification and preliminary SAR studies of (+)-geodin as a glucose uptake stimulator for rat adipocytes. *J. Antibiot. (Tokyo)* **58**(9), 583-589 (2005).

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, 08/23/2022

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