

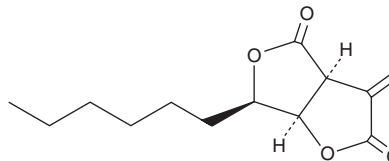
PRODUCT INFORMATION



Sporothriolide

Item No. 29323

CAS Registry No.: 154799-92-5
Formal Name: (3aS,6R,6aR)-6-hexyltetrahydro-3-methylene-furo[3,4-b]furan-2,4-dione
MF: C₁₃H₁₈O₄
FW: 238.3
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Hypoxylon sp.



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

Laboratory Procedures

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

Description

Sporothriolide is a fungal metabolite that has been found in *Sporothrix*.¹ It is active against the plant pathogenic fungi *U. violacea*, *M. microspore*, and *E. repens*, and the alga *C. fusca*, but not the bacteria *B. megaterium* and *E. coli*, when used at a concentration of 10 mg/ml. Sporothriolide inhibits mycelial growth of the plant pathogenic fungus *R. solani* (EC₅₀ = 3.04 µg/ml) and protects against *R. solani*-induced rice sheath blight on rice plants when used at a concentration of 200 µg/ml.² It also prevents necrosis induced by the plant pathogenic fungus *B. cinerea* in pepper seedlings at 500 ppm.¹ Sporothriolide (4 mg/ml) prevents germination of *L. sativum* and *M. sativa* seedlings.

References

1. Krohn, K., Ludewig, K., Aust, H.-J., et al. Biologically active metabolites from fungi. 3. sporothriolide, discosiolide, and 4-*epi*-ethisolide--new furofurandiones from *Sporothrix* sp., *Discosia* sp., and *Pezizula livida*. *J. Antibiot. (Tokyo)* **47**(1), 113-118 (1994).
2. Cao, L.-L., Zhang, Y.-Y., Liu, Y.-J., et al. Anti-phytopathogenic activity of sporothriolide, a metabolite from endophyte *Nodulisporium* sp. A21 in *Ginkgo biloba*. *Pestic. Biochem. Physiol.* **129**, 7-13 (2016).

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, 08/16/2023

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