

PRODUCT INFORMATION



Anhydroophiobolin A

Item No. 25483

CAS Registry No.: 6026-65-9
Formal Name: (2'S,3'S,3aR,5'R,6aS,9aS,10aR)-2,3a,4,4',5',6a,7,9a,10,10a-decahydro-3',9,10a-trimethyl-5'-(2-methyl-1-propen-1-yl)-7-oxo-spiro[dicyclopenta[a,d]cyclooctene-3(1H),2'(3'H)-furan]-6-carboxaldehyde

Synonyms: Anhydrocochliobolin A, 3-Anhydroophiobolin A

MF: C₂₅H₃₄O₃

FW: 382.5

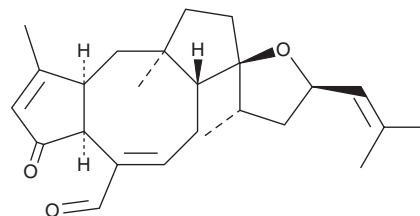
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Bipolaris* sp.



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

Laboratory Procedures

Anhydroophiobolin A is supplied as a solid. A stock solution may be made by dissolving the anhydroophiobolin A in the solvent of choice, which should be purged with an inert gas. Anhydroophiobolin A is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Anhydroophiobolin A is an ophiobolin fungal metabolite that has been found in *C. heterostrophus* fermentation broths.¹ It is cytotoxic to HepG2 and K562 cancer cells (IC₅₀s = 55.7 and 39.5 μM, respectively).²

References

1. Li, E., Clark, A.M., Rotella, D.P., *et al.* Microbial metabolites of ophiobolin A and antimicrobial evaluation of ophiobolins. *J. Nat. Prod.* **58**(1), 74-81 (1995).
2. Wang, Q.-X., Yang, J.-L., Qi, Q.-Y., *et al.* 3-Anhydro-6-hydroxy-ophiobolin A, a new sesterterpene inhibiting the growth of methicillin-resistant *Staphylococcus aureus* and inducing the cell death by apoptosis on K562, from the phytopathogenic fungus *Bipolaris oryzae*. *Bioorg. Med. Chem. Lett.* **23**(12), 3547-3550 (2013).

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, 08/30/2023

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