

PRODUCT INFORMATION



Myceliothermophin E

Item No. 30053

CAS Registry No.: 955083-90-6
Formal Name: (5Z)-1,5-dihydro-5-(2-methylpropylidene)-3-[[[(1S,2R,4aR,6R,8aS)-1,2,4a,5,6,7,8,8a-octahydro-3,4a,6-trimethyl-2-[(1E)-1-methyl-1-propen-1-yl]-1-naphthalenyl]carbonyl]-2H-pyrrol-2-one

MF: C₂₆H₃₇NO₂

FW: 395.6

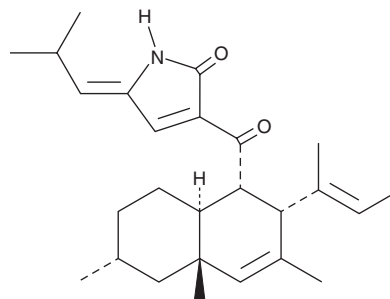
Purity: ≥70%

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years

Item Origin: Fungi/*Myceliophthora* sp.



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

Laboratory Procedures

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

Description

Myceliothermophin E is a polyketide-amino acid hybrid fungal metabolite that has been found in *T. thermophilus* and has anticancer and antimicrobial activities.^{1,2} It is cytotoxic to DLD-1, Hep3B, HepG2, and HGC-27 cancer cells (IC₅₀s = 0.32, 0.42, 0.26, and 0.08 µg/ml, respectively).¹ Myceliothermophin E is active against methicillin-resistant, but not -sensitive, *S. aureus* (MICs = 15.8 and >100 µM, respectively).²

References

1. Gao, Y.-L., Zhang, M.-L., Wang, X., *et al.* Isolation and characterization of a new cytotoxic polyketide-amino acid hybrid from *Thermothelomyces thermophilus* ATCC 42464. *Nat. Prod. Res.* **Epub ahead of print**, (2019).
2. Wang, X., Zhao, L., Liu, C., *et al.* New tetramic acids comprising of decalin and pyridones from *Chaetomium olivaceum* SD-80A with antimicrobial activity. *Front. Microbiol.* **10**, 2958 (2020).

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, 05/20/2020

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