

PRODUCT INFORMATION



Cyclothiocurvularin A

Item No. 31668

Formal Name: (1R,3aR,7S,15aR)-1,12,14-trihydroxy-7-methyl-9,15-dioxo-1,2,3a,4,6,7,9,10,15,15a-decahydro-5H-benzo[d]thieno[3,2-g][1]oxacyclododecine-1-carboxylic acid

MF: C₁₉H₂₂O₈S

FW: 410.4

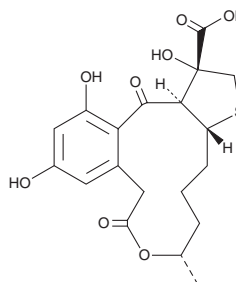
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Curvularia* sp.



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

Laboratory Procedures

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

Description

Cyclothiocurvularin A is a macrocyclic polyketide that has been found in *Penicillium*.¹

Reference

1. de Castro, M.V., Ióca, L.P., Williams, D.E., *et al.* Condensation of macrocyclic polyketides produced by *Penicillium* sp. DRF2 with mercaptopyruvate represents a new fungal detoxification pathway. *J. Nat. Prod.* **79**(6), 1668-1678 (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, 08/11/2023

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