

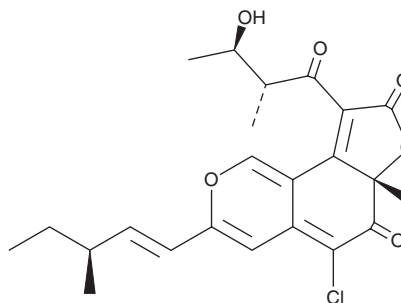
PRODUCT INFORMATION



Chaetoviridin A

Item No. 31174

CAS Registry No.: 128252-98-2
Formal Name: (6aS)-5-chloro-9-[(2R,3R)-3-hydroxy-2-methyl-1-oxobutyl]-6a-methyl-3-[[1E,3S)-3-methyl-1-penten-1-yl]-6H-furo[2,3-h]-2-benzopyran-6,8(6aH)-dione
MF: C₂₃H₂₅ClO₆
FW: 432.9
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Chaetomium* sp.



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

Laboratory Procedures

Chaetoviridin A is supplied as a solid. A stock solution may be made by dissolving the chaetoviridin A in the solvent of choice, which should be purged with an inert gas. Chaetoviridin A is soluble in organic solvents such as acetone, chloroform, DMSO, and methanol.

Description

Chaetoviridin A is a fungal metabolite that has been found in *Chaetomium*.¹ Plant pathogenic, antitumor, and cholesteryl ester transfer protein (CETP) inhibitory activity attributed to chaetoviridin A was determined prior to its structural reassignment by total synthesis, NMR, and circular dichroism analysis.¹⁻⁴

References

1. Makrourgras, M., Coffinier, R., Oger, S., *et al.* Total synthesis and structural revision of chaetoviridins A. *Org. Lett.* **19(15)**, 4146-4149 (2017).
2. Park, J.-H., Choi, G.J., Jang, K.S., *et al.* Antifungal activity against plant pathogenic fungi of chaetoviridins isolated from *Chaetomium globosum*. *FEMS Microbiol. Lett.* **252(2)**, 309-313 (2005).
3. Yasukawa, K., Takahashi, M., Natori, S., *et al.* Azaphilones inhibit tumor promotion by 12-O-tetradecanoylphorbol-13-acetate in two-stage carcinogenesis in mice. *Oncology* **51(1)**, 108-112 (1994).
4. Tomoda, H., Matsushima, C., Tabata, N., *et al.* Structure-specific inhibition of cholesteryl ester transfer protein by azaphilones. *J. Antibiot. (Tokyo)* **52(2)**, 160-170 (1999).

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, 08/10/2023

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