

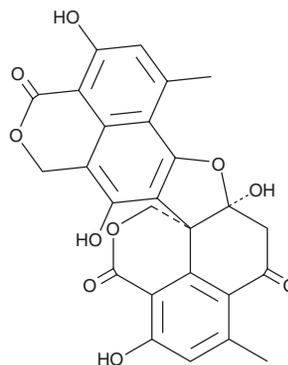
PRODUCT INFORMATION



Bacillosporin C

Item No. 23877

CAS Registry No.: 76706-63-3
Formal Name: 5,5a-dihydro-1,5a,9,13-tetrahydro-3,7-dimethyl-4H,10H,12H,14H,16H-dibenzo[de,d'e']furo[2,3-g:5,4-i']bis[2]benzopyran-4,10,16-trione
MF: C₂₆H₁₈O₁₀
FW: 490.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: *Aspergillus* sp.



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

Laboratory Procedures

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

Description

Bacillosporin C is an oxaphenalenone dimer originally isolated from *T. bacillosporus*.¹ Bacillosporin C, an anhydride, is formed from the lactone bacillosporin D in the mangrove endophytic fungus SBE-14.² Similar oxaphenalenone dimers have antibiotic activity and inhibit acetylcholinesterase.³

References

1. Yamazaki, M. and Okuyama, E. Isolation and structures of oxaphenalenone dimers from *Talaromyces bacillosporus*. *Chem. Pharm. Bull.* **28(12)**, 3649-3655 (1980).
2. Guo, Z., Shao, C., She, Z., et al. ¹H and ¹³C NMR assignments for two oxaphenalenones bacillosporin C and D from the mangrove endophytic fungus SBE-14. *Magn. Reson. Chem.* **45(5)**, 439-441 (2007).
3. Wu, B., Ohlendorf, B., Oesker, V., et al. Acetylcholinesterase inhibitors from a marine fungus *Talaromyces* sp. strain LF458. *Mar. Biotechnol. (NY)* **17(1)**, 110-119 (2015).

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, 11/16/2022

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