

PRODUCT INFORMATION



Sesquicillin A

Item No. 29326

CAS Registry No.: 51103-58-3
Formal Name: 3-[[[(1S,4aS,5R,6R,8aS)-6-(acetyloxy)decahydro-5,8a-dimethyl-2-methylene-5-(4-methyl-3-penten-1-yl)-1-naphthalenyl]methyl]-4-hydroxy-5,6-dimethyl-2H-pyran-2-one

Synonym: Sesquicillin

MF: C₂₉H₄₂O₅

FW: 470.6

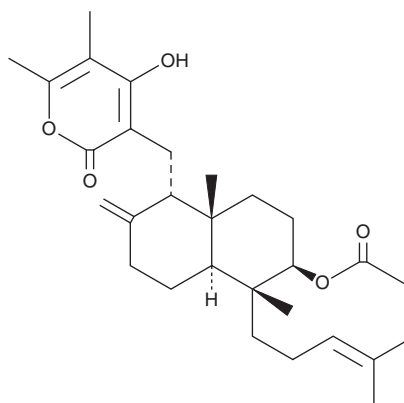
Purity: ≥70%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

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

Description

Sesquicillin A is a fungal metabolite that has been found in *Albophoma*.¹ It inhibits the growth of *A. salina* brine shrimp (MIC = 6.25 µg/ml). Sesquicillin is cytotoxic to Jurkat cells (IC₅₀ = 34 µM). It induces cell cycle arrest at the G₁ phase and inhibits proliferation of MCF-7 breast cancer cells when used at a concentration of 20 µg/ml.²

References

1. Uchida, R., Imasato, R., Yamaguchi, Y., *et al.* New sesquicillins, insecticidal antibiotics produced by *Albophoma sp.* FKI-1778. *J. Antibiot. (Tokyo)* **58(6)**, 397-404 (2005).
2. Jeong, H.W., Lee, H.J., Kho, Y.H., *et al.* Biological effects of G₁ phase arrest compound, sesquicillin, in human breast cancer cell lines. *Bioorg. Med. Chem.* **10(10)**, 3129-3134 (2002).

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, 08/16/2023

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