

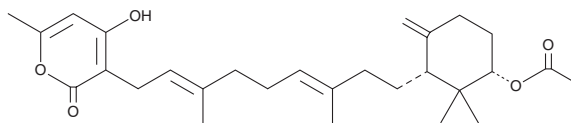
PRODUCT INFORMATION



Sartorypyrone A

Item No. 23863

CAS Registry No.: 1452396-10-9
Formal Name: *rel-(+)-3-[(2E,6E)-9-[(1R,3S)-3-(acetyloxy)-2,2-dimethyl-6-methylenecyclohexyl]-3,7-dimethyl-2,6-nonadien-1-yl]-4-hydroxy-6-methyl-2H-pyran-2-one*
Synonym: 20-O-Acetyl Sartorypyrone D
MF: C₂₈H₄₀O₅
FW: 456.6
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Neosartorya* sp.



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

Laboratory Procedures

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

Description

Sartorypyrone A is a meroditerpene metabolite produced by *Neosartorya* fungal species.¹ It inhibits the growth of MCF-7 breast, NCI-H460 lung, and A375-C5 melanoma cancer cell lines (GI₅₀s = 46.3, 37.3, and 21.5 μM, respectively). Sartorypyrone A inhibits growth of Gram-positive bacteria, including multidrug-resistant *S. aureus* (MIC = 32 μg/ml), and reduces the mass of *S. aureus* and *B. subtilis* biofilms *in vitro*.²

References

1. Eamvijarn, A., Gomes, N.M., Dethoup, T., *et al.* Bioactive meroditerpenes and indole alkaloids from the soil fungus *Neosartorya fischeri* (KUFC 6344), and the marine-derived fungi *Neosartorya lacinosia* (KUFC 7896) and *Neosartorya tsunodae* (KUFC 9213). *Tetrahedron* **69(40)**, 8583-8591 (2013).
2. Gomes, N.M., Bessa, L.J., Buttachon, S., *et al.* Antibacterial and antibiofilm activities of tryptoquivalines and meroditerpenes isolated from the marine-derived fungi *Neosartorya paulistensis*, *N. lacinosia*, *N. tsunodae*, and the soil fungi *N. fischeri* and *N. siamensis*. *Mar. Drugs* **12(2)**, 822-839 (2014).

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, 08/23/2023

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