

PRODUCT INFORMATION



Podosporin A

Item No. 44804

CAS Registry No.: 176018-99-8

Formal Name: [3 α ,4 α β ,7 α ,7 α ,11(1R*,3E),13 α S*]-
1,2,3,4,4 α ,5,6,7,7 α ,8-decahydro-
3-hydroxy-4,4,7,7 α -tetramethyl-
10-(1-methyl-2-oxo-3-pentenyl)-
benzo[d]xanthene-9,12-dione
($\pm$)-Podosporin A

Synonym:

MF: C₂₇H₃₆O₅

FW: 440.6

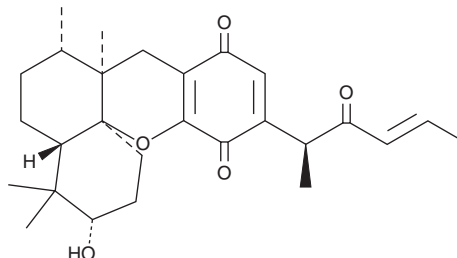
Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Fungus/*Podospira* sp.



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

Laboratory Procedures

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

Description

Podospirin A is a sesquiterpene quinone and fungal metabolite that has been found in *P. decipiens* and has diverse biological activities.¹ It is active against *S. fimicola* and *A. furfuraceus* (MICs = 175 and 50 $\mu\text{g/ml}$, respectively), as well as *B. subtilis*, *S. aureus*, and *C. albicans* when used in a disc assay at a concentration of 50 $\mu\text{g/disc}$. Podospirin A is toxic to brine shrimp (LC_{50} = 10 $\mu\text{g/ml}$).

Reference

- 1 Weber, H.A., Baenziger, N.C., and Gloer, J.B. Podospirin A: A novel antifungal metabolite from the coprophilous fungus *Podospira decipiens* (Wint.) niessl. *J. Org. Chem.* **53**(19), 4567-4569 (1988).

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, 11/24/2025

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