

PRODUCT INFORMATION



Aranorosin

Item No. 27478

CAS Registry No.: 117184-53-9
Formal Name: (2E,4E,6R)-N-[(1 α ,2 β ,3 α ,4' β ,5 α ,5'R,7 α)-dihydro-5'-hydroxy-6-oxospiro[4,8-dioxatricyclo[5.1.0.0^{3,5}]octane-2,2'(3'H)-furan]-4'-yl]-4,6-dimethyl-2,4-dodecadienamide

MF: C₂₃H₃₃NO₆

FW: 419.5

Purity: $\geq 95\%$

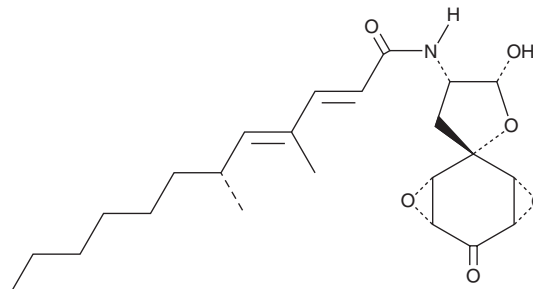
UV/Vis.: λ_{\max} : 265 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Fungus/*Gymnascella* sp.



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

Laboratory Procedures

Aranorosin is supplied as a solid. A stock solution may be made by dissolving the aranorosin in the solvent of choice, which should be purged with an inert gas. Aranorosin is soluble in the organic solvent ethanol at a concentration of approximately 1 mg/ml. Aranorosin is soluble in methanol and DMSO.

Description

Aranorosin is a fungal metabolite originally isolated from *P. roseus*.¹ It has antimicrobial activity against *B. subtilis*, *A. niger*, and *C. albicans* when used at a concentration of 1 mg/ml. Aranorosin also reduces viability in apoptosis-resistant HeLa/Bcl-2 cells.²

References

1. Vijayakumar, E.K.S., Roy, K., Chatterjee, S., *et al.* Structure-activity relationships of new aranorosin analogs to antifungal activity. *J. Antibiot. (Tokyo)* **51(5)**, 522-524 (1998).
2. Nakashima, T., Tanaka, R., Yamashita, Y., *et al.* Aranorosin and a novel derivative inhibit the anti-apoptotic functions regulated by Bcl-2. *Biochem. Biophys. Res. Commun.* **377(4)**, 1085-1090 (2008).

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, 08/09/2023

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