

PRODUCT INFORMATION



Agonodepside B

Item No. 32930

CAS Registry No.: 445298-41-9
Formal Name: 2,4-dihydroxy-3-methyl-6-[(1E)-1-methyl-1-propen-1-yl]-benzoic acid, 4-carboxy-3-hydroxy-2-methyl-5-[(1E)-1-methyl-1-propen-1-yl]phenyl ester

MF: C₂₄H₂₆O₇

FW: 426.5

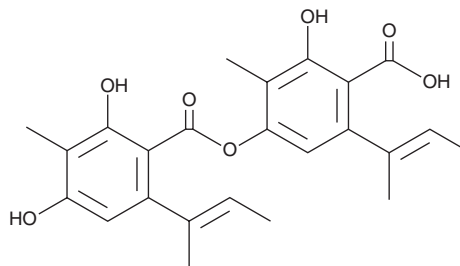
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungi/*Emericella* sp.



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

Laboratory Procedures

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

Description

Agonodepside B is a fungal metabolite originally isolated from the endophytic fungus F7524.¹

Reference

1. Cao, S., Lee, A.S.Y., Huang, Y., *et al.* Agonodepsides A and B: Two new depsides from a filamentous fungus F7524. *J. Nat. Prod.* **65**(7), 1037-1038 (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, 11/10/2022

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