

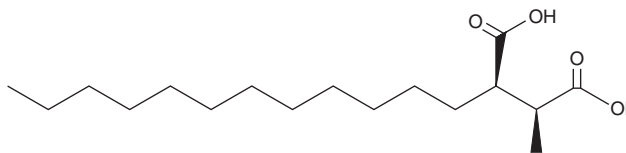
PRODUCT INFORMATION



Roccellic Acid

Item No. 29909

CAS Registry No.: 29838-46-8
Formal Name: 2R-dodecyl-3S-methylbutanedioic acid
Synonyms: (+)-Roccellic Acid, NSC 249985
MF: C₁₇H₃₂O₄
FW: 300.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Lecanora* sp.



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

Laboratory Procedures

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

Description

Roccellic acid is a lichen secondary metabolite that has been found in *R. montagnei* and has antibacterial and anticancer activities.^{1,2} It is active against the bacteria *S. gordonii* and *P. gingivalis* (MIC = 46.9 µg/ml for both).¹ Roccellic acid (100 µg/ml) inhibits proliferation of MDA-MB-231, MCF-7, and DLD-1 cancer cells by 65.3, 75.8, and 87.9%, respectively.²

References

1. Sweidan, A., Chollet-Krugler, M., Sauvager, A., *et al.* Antibacterial activities of natural lichen compounds against *Streptococcus gordonii* and *Porphyromonas gingivalis*. *Fitoterapia* **121**, 164-169 (2017).
2. Mishra, T., Shukla, S., Meena, S., *et al.* Isolation and identification of cytotoxic compounds from a fruticose lichen *Roccella montagnei*, and it's *in silico* docking study against CDK-10. *Rev. Bras. Farmacogn.* **27(6)**, 724-728 (2017).

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