

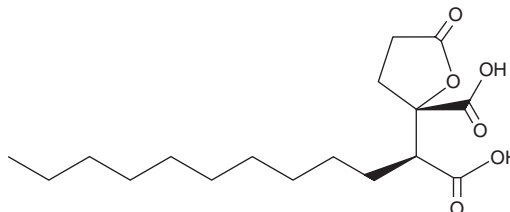
PRODUCT INFORMATION



Spiculisporic Acid

Item No. 31334

CAS Registry No.: 469-77-2
Formal Name: (αS,2S)-2-carboxy-α-decyltetrahydro-5-oxo-2-furanacetic acid
Synonyms: (-)-Spiculisporic Acid, NSC 244387
MF: C₁₇H₂₈O₆
FW: 328.4
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Spiculisporic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the spiculisporic acid in the solvent of choice, which should be purged with an inert gas. Spiculisporic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of spiculisporic acid in these solvents is approximately 30 mg/ml.

Description

Spiculisporic acid is a biosurfactant that has been found in *P. spiculisporum*.¹

Reference

1. Ishigami, Y., Zhang, Y. and Goto, M. Molecular and crystal structure of spiculisporic acid and correlation with the surface activity. *J. Oleo Sci.* **62**(10), 795-802 (2013).

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, 12/09/2022

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