

PRODUCT INFORMATION



Berkeleylactone E

Item No. 29826

CAS Registry No.: 122211-62-5
Formal Name: [5R-(3E,5R,6S,16R)]-butanedioic acid, mono(5-hydroxy-16-methyl-2-oxooxacyclohexadec-3-en-6-yl) ester

MF: C₂₀H₃₂O₇

FW: 384.5

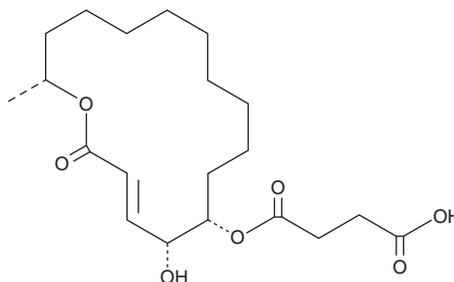
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Penicillium* sp.



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

Laboratory Procedures

Berkeleylactone E is supplied as a solid. A stock solution may be made by dissolving the berkeleylactone E in the solvent of choice, which should be purged with an inert gas. Berkeleylactone E is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

Berkeleylactone E is a macrolide antibiotic that has been found in *P. fuscum* and *P. camembertii/clavigerum* co-culture.¹ It is active against *S. aureus* (MIC = 125 μM).

Reference

1. Stierle, A.A., Stierele, D.B., Decato, D., *et al.* The berkeleylactones, antibiotic macrolides from fungal coculture. *J. Nat. Prod.* **80**(4), 1150-1160 (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, 08/10/2023

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