

PRODUCT INFORMATION



Malolactomycin C

Item No. 31673

CAS Registry No.: 189759-03-3
Formal Name: (+)-1-[5,7,9,11,21,25,27,29,31,35,37,38,39-tridecahydroxy-17-[9-[[imino(methylamino)methyl]amino]-1,3-dimethyl-3-nonen-1-yl]-10,16,20,22,26,30,34-heptamethyl-19-oxo-18,41-dioxabicyclo[35.3.1]hentetraconta-12,14,22-trien-3-yl] ester, propanedioic acid

MF: C₆₂H₁₀₉N₃O₂₀

FW: 1,216.6

Purity: ≥95%

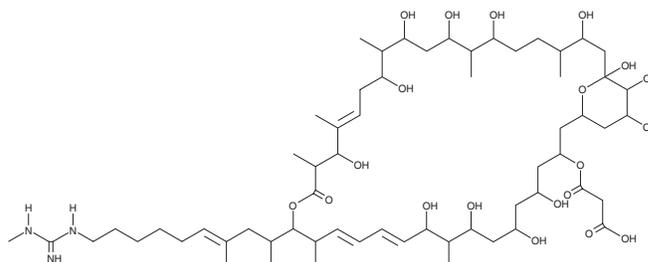
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Streptomyces* sp.

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



Laboratory Procedures

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

Description

Malolactomycin C is a bacterial metabolite originally isolated from *Streptomyces* that has fungicidal activity.¹ It is active against the plant pathogenic fungi *B. cinerea*, *C. fulvum*, and *P. oryzae* (MIC = 25 µg/ml for all).

Reference

1. Tanaka, Y., Yoshida, H., Enomoto, Y., *et al.* Malolactomycins C and D, new 40-membered macrolides active against *Botrytis*. *J. Antibiot. (Tokyo)* **50(3)**, 194-200 (1997).

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