

PRODUCT INFORMATION



Martinomycin

Item No. 27737

CAS Registry No.: 160791-16-2

Formal Name: (α R,2S,3S,4S,5S,6S)-6-[[[(2R,3R,4R,5S,7S,9R,10R)-3,9-dimethoxy-2,4,10-trimethyl-2-[(2S,2'R,5R,5'R)-octahydro-5'-[(2S,3S,4S,5R,6S)-tetrahydro-6-hydroxy-3,5,6-trimethyl-4-[[[(2R,5S,6R)-tetrahydro-5-methoxy-6-methyl-2H-pyran-2-yl]oxy]-2H-pyran-2-yl][2,2'-bifuran]-5-yl]-1,6-dioxaspiro[4.5]dec-7-yl]methyl]tetrahydro-2-hydroxy-4,5-dimethoxy- α ,3,5-trimethyl-2H-pyran-2-acetic acid

Synonym: LL-D 37187 α

MF: C₄₉H₈₄O₁₇

FW: 945.2

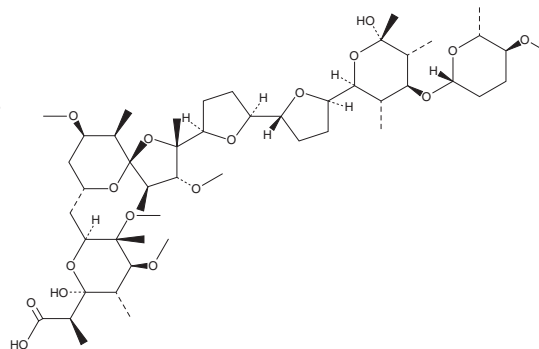
Purity: \geq 95%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Bacterium/*Actinoallamurus yoronensis*



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

Laboratory Procedures

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

Description

Martinomycin is a polyether antibiotic originally isolated from the actinomycete *S. salivialis*.^{1,2} It is active against a variety of bacteria, including 11 strains of *S. aureus*, three strains of *S. pneumoniae*, and *E. faecalis* (MICs = 0.12-0.5, 0.06, and 0.5 μ g/ml, respectively).²

References

1. Carter, G.T., Schlingmann, G., Kenion, G.B., *et al.* Martinomycin, a new polyether antibiotic produced by *Streptomyces salivialis*. II. Isolation and structure determination. *J. Antibiot. (Tokyo)* **47(12)**, 1549-1553 (1994).
2. Bernan, V.S., Montenegro, D.A., Goodman, J.J., *et al.* Martinomycin, a new polyether antibiotic produced by *Streptomyces salivialis*. I. Taxonomy, fermentation and biological activity. *J. Antibiot. (Tokyo)* **47(12)**, 1434-1441 (1994).

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, 08/25/2023

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