

PRODUCT INFORMATION



Reveromycin A

Item No. 17458

CAS Registry No.: 134615-37-5
Formal Name: butanedioic acid, 1-[(2S,3R,6S,8R,9S)-3-butyl-8-[(2E,4E,6S,7S,8E)-9-carboxy-6-hydroxy-3,7-dimethyl-2,4,8-nonatrien-1-yl]-2-[(1E,3E)-4-carboxy-3-methyl-1,3-butadien-1-yl]-9-methyl-1,7-dioxaspiro[5.5]undec-3-yl] ester

MF: C₃₆H₅₂O₁₁

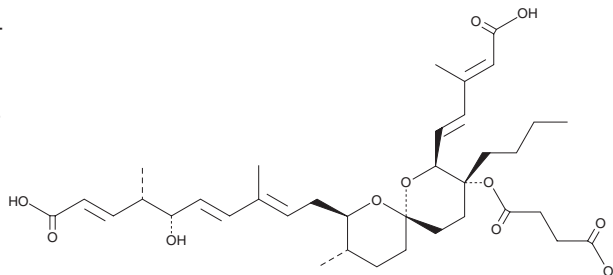
FW: 660.8

Purity: ≥99%

Supplied as: A lyophilisate

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

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

Description

Reveromycin A is the major component of a complex of spiroketal antibiotics isolated from *Streptomyces* sp. It inhibits the mitogenic activity of epidermal growth factor in Balb/MK cells (IC₅₀ = 0.7 µg/ml), displays antiproliferative activity against human KB and K562 tumor cell lines (IC₅₀s = 1.9 and 1.6 µg/ml, respectively), and demonstrates antifungal activity against *C. albicans* (MIC = 2 µg/ml at pH 3).¹ Reveromycin A also has been shown to inhibit bone resorption by inducing apoptosis in osteoclasts with an IC₅₀ value of 0.7 µM.²

References

1. Takahashi, H., Osada, H., Koshino, H., *et al.* Reveromycins, new inhibitors of eukaryotic cell growth. II. Biological activities. *J. Antibiot. (Tokyo)* **45(9)**, 1414-1419 (1992).
2. Woo, J.-T., Kawatani, M., Kato, M., *et al.* Reveromycin A, an agent for osteoporosis, inhibits bone resorption by inducing apoptosis specifically in osteoclasts. *Proc. Natl. Acad. Sci. USA* **103(12)**, 4729-4734 (2006).

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, 12/12/2022

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