

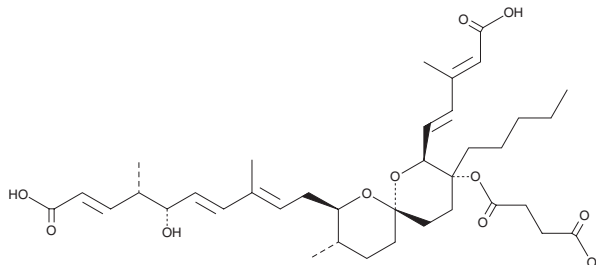
# PRODUCT INFORMATION



## Reveromycin D

Item No. 28101

**CAS Registry No.:** 144860-70-8  
**Formal Name:** butanedioic acid, 1-[(2S,3R,6S,8R,9S)-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-3-pentyl-1,7-dioxaspiro[5.5]undec-3-yl] ester  
**MF:** C<sub>37</sub>H<sub>54</sub>O<sub>11</sub>  
**FW:** 674.8  
**Purity:** ≥95%  
**Supplied as:** A lyophilisate  
**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

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

Reveromycin D is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Reveromycin D is a bacterial metabolite originally isolated from *Streptomyces*.<sup>1</sup> It inhibits EGF-induced mitogenic activity in Balb/MK cells and has pH-dependent antifungal activity against *C. albicans* (MICs = 2 and >500 µg/ml at pH 3 and 7.4, respectively).<sup>2</sup> Reveromycin D also inhibits proliferation of KB and K562 cells (IC<sub>50</sub>s = 1.6 and 1.3 µg/ml, respectively).

### References

1. Takahashi, H., Osada, H., Koshino, H., *et al.* Reveromycins, new inhibitors of eukaryotic cell growth. I. Producing organism, fermentation, isolation and physico-chemical properties. *J. Antibiot. (Tokyo)* **45(9)**, 1409-1413 (1992).
2. 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).

#### 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, 10/11/2022

#### 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