

PRODUCT INFORMATION

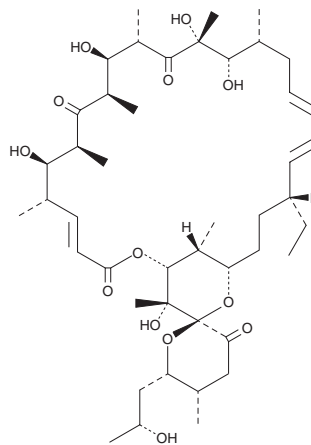


Oligomycin E

Item No. 20185

CAS Registry No.: 110231-34-0
Formal Name: (1R,2'S,4E,5'S,6S,6'S,7R,8S,10R,11R,12S,14R,15S,16R,18E,20E,22R,25S,28R,29R)-22-ethyl-5',6'-dihydro-7,11,14,15,28-pentahydroxy-6'-[(2R)-2-hydroxypropyl]-5',6,8,10,12,14,16,28,29-nonamethyl-spiro[2,26-dioxabicyclo[23.3.1]nonacos-4,18,20-triene-27,2'-[2H]pyran]-3,3',9,13(4'H)-tetrone
26-hydroxy-28-oxo-Oligomycin A

Synonym:
MF: C₄₅H₇₂O₁₃
FW: 821.1
Purity: ≥95%
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

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

Oligomycin E 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

Oligomycin E is a minor metabolite of the oligomycin complex (Item No. 11341) produced by several species of *Streptomyces*. It exhibits relatively weak antifungal activity compared to other oligomycins, yet is active against Gram-positive bacteria and demonstrates strong antitumor activity against HeLa cells (IC₅₀ = 14 ng/ml).¹

Reference

1. Kobayashi, K., Nishino, C., Ohya, J. *et al.* Oligomycin E, a new antitumor antibiotic produced by *Streptomyces* sp. MCI-2225. *J. Antibiot. (Tokyo)* **40(7)**, 1053-1057 (1987).

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, 12/15/2022

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