

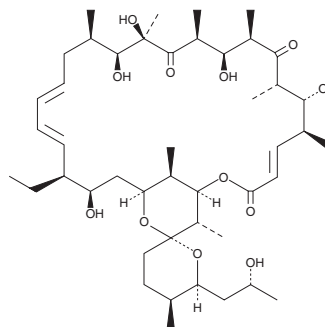
PRODUCT INFORMATION



21-hydroxy Oligomycin A

Item No. 20599

CAS Registry No.: 102042-09-1
Formal Name: (1R,2'S,4E,5'S,6S,6'S,7R,8S,10R,11R,12S,14R,15S,16R,18E,20E,22R,25S,28S,29R)-22-ethyl-3',4',5',6'-tetrahydro-7,11,14,15,23-pentahydroxy-6'-[(2R)-2-hydroxypropyl]-5',6,8,10,12,14,16,28,29-nonamethyl-spiro[2.26-dioxabicyclo[23.3.1]nonacosa-4,18,20-triene-27,2'-[2H]pyran]-3,9,13-trione



MF: C₄₅H₇₄O₁₂

FW: 807.1

Purity: ≥95%

Supplied as: A lyophilisate

Storage: -20°C

Stability: ≥2 years

Item Origin: Bacterium/*Streptomyces* sp.

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

Laboratory Procedures

21-hydroxy Oligomycin A is supplied as a lyophilisate. A stock solution may be made by dissolving the 21-hydroxy oligomycin A in the solvent of choice, which should be purged with an inert gas. 21-hydroxy Oligomycin A is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

21-hydroxy Oligomycin A is an antibiotic that was first isolated from *S. cyaneogriseus* ssp. *noncyanogenus* (LL-F28249). It is reported to be cytotoxic to human colon cancer SW620 cells (IC₅₀ = 14.4 μM), to inhibit the ABC transporter efflux pump P-glycoprotein, and to prevent K-Ras plasma membrane localization (IC₅₀ = 4.82 nM).¹

Reference

1. Salim, A.A., Tan, L., Huang, X.-C., *et al.* Oligomycins as inhibitors of K-Ras plasma membrane localisation. *Org. Biomol. Chem.* **14**(2), 711-715 (2016).

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, 01/15/2021

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