

PRODUCT INFORMATION



Heliquinomycin

Item No. 32922

CAS Registry No.: 178182-49-5
Formal Name: (2R,3R,3'R)-3'-[(2,6-dideoxy-3-O-methyl- α -L-ribo-hexopyranosyl)oxy]-4,5',8',9-tetrahydro-3,4',9',10-tetrahydroxy-7'-methoxy-5',8',9-trioxo-spiro[benzo[1,2-b:5,4-c']dipyran-2(3H),2'(3'H)-naphtho[2,3-b]furan]-7-carboxylic acid, methyl ester

Synonyms: NSC 702208, Rubymycin

MF: C₃₃H₃₀O₁₇

FW: 698.6

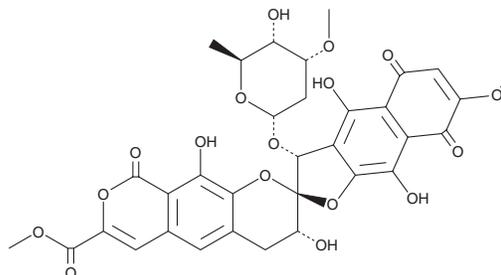
Purity: \geq 80%

Supplied as: A solid

Storage: 4°C

Stability: \geq 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

Heliquinomycin is supplied as a solid. A stock solution may be made by dissolving the heliquinomycin in the solvent of choice, which should be purged with an inert gas. Heliquinomycin is soluble in ethanol and DMSO.

Description

Heliquinomycin is a bacterial metabolite originally isolated from *Streptomyces* that has diverse biological activities.¹ It is active against a variety of Gram-positive bacteria, including strains of *B. anthracis*, *B. subtilis*, and methicillin-sensitive or -resistant *S. aureus* (MICs = <0.05-0.39 μ g/ml). Heliquinomycin inhibits the activity of DNA helicase with a K_i value of 6.8 μ M. It reduces the growth of L1210 leukemia, B16 melanoma, and FS-3 fibrosarcoma cells (IC₅₀s = 0.97, 0.89, and 0.83 μ g/ml, respectively).

Reference

1. Chino, M., Nishikawa, K., Umekita, M., *et al.* Heliquinomycin, a new inhibitor of DNA helicase, produced by *Streptomyces* sp. MJ929-SF2 I. Taxonomy, production, isolation, physico-chemical properties and biological activities. *J. Antibiot. (Tokyo)* **49(8)**, 752-757 (1996).

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, 03/07/2023

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