

PRODUCT INFORMATION



4'-Acetyl Chrysomycin B

Item No. 33024

Formal Name: (2R,3S,4S,5S,6S)-4,5-dihydroxy-6-(1-hydroxy-10,12-dimethoxy-6-oxo-8-vinyl-6H-dibenzo[c,h]chromen-4-yl)-2,4-dimethyltetrahydro-2H-pyran-3-yl acetate

MF: C₃₀H₃₀O₁₀

FW: 550.6

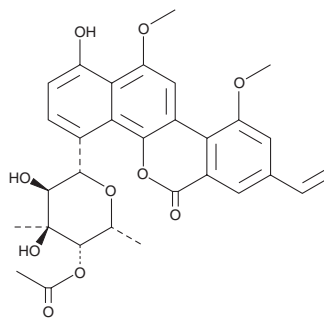
Purity: ≥95%

Supplied as: A solid

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

4'-Acetyl chrysomycin B is supplied as a solid. A stock solution may be made by dissolving the 4'-acetyl chrysomycin B in the solvent of choice, which should be purged with an inert gas. 4'-Acetyl chrysomycin B is soluble in methanol and DMSO.

Description

4'-Acetyl chrysomycin B is a bacterial metabolite and derivative of chrysomycin B (Item No. 14189) that has been found in *Streptomyces* and has antibacterial and anticancer activities.¹ It is active against strains of *B. subtilis*, *B. cereus*, *Corynebacterium bovis*, and *E. faecalis* (MICs = 2-64 µg/ml). 4'-Acetyl chrysomycin B is cytotoxic against a panel of human cancer cell lines, including doxorubicin-sensitive or -resistant cells (IC₅₀s = 1.7-14 and 260-370 ng/ml, respectively).

Reference

1. Wada, S.-I., Sawa, R., Iwanami, F., *et al.* Structures and biological activities of novel 4'-acetylated analogs of chrysomycins A and B. *J. Antibiot. (Tokyo)* **70(11)**, 1078-1082 (2017).

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, 07/28/2023

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