

PRODUCT INFORMATION



Saccharocarcin A

Item No. 28094

CAS Registry No.: 158475-32-2
Formal Name: 10-[[4-(acetylamino)-2,4,6-trideoxy-3-C-methylhexopyranosyl]oxy]-4-[[O-2,6-dideoxyhexopyranosyl-(1→4)-O-2,6-dideoxy-3-O-(tetrahydro-5-hydroxy-6-methyl-2H-pyran-2-yl)hexopyranosyl-(1→4)-2,6-dideoxyhexopyranosyl]oxy]-15-ethyl-2,3,4,4a,6a,9,10,12a,15,16,20a,20b-dodecahydro-21-hydroxy-1,3,7,9,11,20a-hexamethyl-18H-16a,19-metheno-16aH-benzo[b]naphth[2,1-j]oxacyclotetradecin-18,20(1H)-dione

MF: C₆₇H₁₀₁NO₂₀
FW: 1,240.5

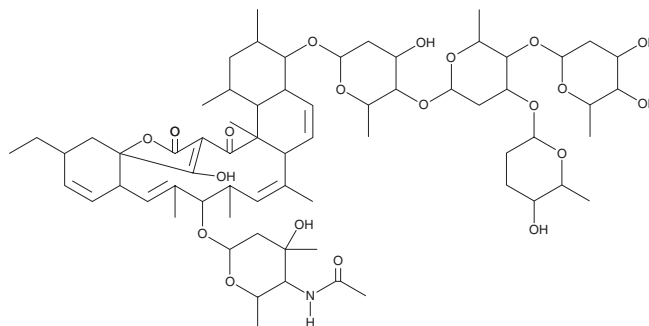
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Amycolatopsis* sp.



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

Laboratory Procedures

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

Description

Saccharocarcin A is an antibiotic originally isolated from *S. aerocolonigenes* subsp. *antibiotica*.¹ It is active against *M. luteus*, *S. aureus*, *E. coli*, *P. aeruginosa*, and *C. albicans* in a disc assay. Saccharocarcin A inhibits *C. trachomatis* infection by 88% without inducing cytotoxicity in McCoy cells when used at a concentration of 0.5 µg/ml.

Reference

1. Horan, A.C., Shearer, M.C., Hegde, V., et al. A family of novel macrocyclic lactones, the saccharocarmins produced by *Saccharothrix aerocolonigenes* subsp. *antibiotica*. I. Taxonomy, fermentation, isolation and biological properties. *J. Antibiot. (Tokyo)* **50(2)**, 119-125 (1997).

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, 11/01/2022

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