

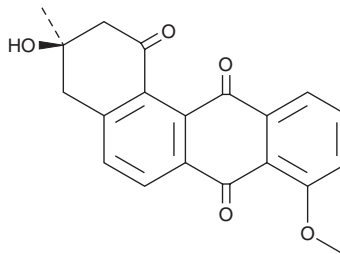
PRODUCT INFORMATION



6-Deoxy-8-O-methylrabelomycin

Item No. 40662

CAS Registry No.: 117620-87-8
Formal Name: 3R,4-dihydro-3-hydroxy-8-methoxy-3-methyl-benz[a]anthracene-1,7,12(2H)-trione
Synonyms: (-)-8-O-Methyltetrangomycin, MM 47755
MF: C₂₀H₁₆O₅
FW: 336.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp. Gö 40/14



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

Laboratory Procedures

6-Deoxy-8-O-methylrabelomycin is supplied as a solid. A stock solution may be made by dissolving the 6-deoxy-8-O-methylrabelomycin in the solvent of choice, which should be purged with an inert gas. 6-Deoxy-8-O-methylrabelomycin is slightly soluble (0.1-1 mg/ml) in acetonitrile and methanol.

Description

6-Deoxy-8-O-methylrabelomycin is an anthraquinone that has been found in *Pseudonocardia* with antiparasitic and antibiotic activities.^{1,2} It is active against the parasite *P. berghei* (IC₅₀ = 18.5 μM) and the bacterium *B. subtilis* (MIC = 25 μg/ml) but does not induce cellular toxicity in HepG2 cells (IC₅₀ = >100 μM).¹ 6-Deoxy-8-O-methylrabelomycin is also active against methicillin-resistant *S. aureus* (MRSA; IC₅₀ = 48 μg/ml) and reduces MRSA production of the virulence factor staphyloxanthin (IC₅₀ = 6 μg/ml).²

References

1. Carr, G., Derbyshire, E.R., Caldera, E., *et al.* Antibiotic and antimalarial quinones from fungus-growing ant-associated *Pseudonocardia* sp. *J. Nat. Prod.* **75(10)**, 1806-1809 (2012).
2. Sakai, K., Koyama, N., Fukuda, T., *et al.* Search method for inhibitors of staphyloxanthin production by methicillin-resistant *Staphylococcus aureus*. *Biol. Pharm. Bull.* **35(1)**, 48-53 (2012).

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, 04/22/2024

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