

PRODUCT INFORMATION

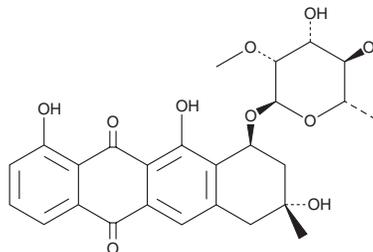


Aranciamycin A

Item No. 31659

CAS Registry No.: 960622-70-2
Formal Name: (8R,10S)-10-[[6-deoxy-2-O-methyl- α -L-mannopyranosyl]oxy]-7,8,9,10-tetrahydro-1,8,11-trihydroxy-8-methyl-5,12-naphthacenedione

MF: C₂₆H₂₈O₁₀
FW: 500.5
Purity: \geq 95%
UV/Vis.: λ_{max} : 226, 260, 433 nm
Supplied as: A solid
Storage: -20°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

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

Description

Aranciamycin A is an anthracyclic antibiotic that has been found in *Streptomyces*.¹ It is active against *M. bovis* and *B. subtilis* (MICs = 30 and 7.5 μ M, respectively) but not *S. aureus*, *E. coli*, *P. aeruginosa*, or *C. albicans* (MICs = $>$ 30 μ M for all).

Reference

1. Khalil, Z.G., Raju, R., Piggott, A.M., et al. Aranciamycins I and J, antimycobacterial anthracyclines from an Australian marine-derived *Streptomyces* sp. *J. Nat. Prod.* **78(4)**, 949-952 (2015).

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, 12/02/2022

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