

# PRODUCT INFORMATION



## Aranciamycin

Item No. 29575

**CAS Registry No.:** 72389-06-1  
**Formal Name:** (2S,3S,4R)-4-[(6-deoxy-2-O-methyl- $\alpha$ -L-mannopyranosyl)oxy]-3,4-dihydro-2,5,7-trihydroxy-3-methoxy-2-methyl-1,6,11(2H)-naphthacenetriene

**Synonym:** NSC 369226

**MF:** C<sub>27</sub>H<sub>28</sub>O<sub>12</sub>

**FW:** 544.5

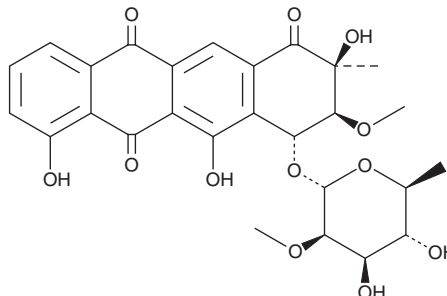
**Purity:**  $\geq$ 70%

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

### Description

Aranciamycin is a fungal metabolite that has been found in *Streptomyces* and has diverse biological activities.<sup>1,2</sup> It inhibits *C. histolyticum* collagenase but not elastase or trypsin in cell-free assays (IC<sub>50</sub>s = 0.37, >10, and >10  $\mu$ M, respectively).<sup>1</sup> Aranciamycin is active against *M. bovis* and *B. subtilis* (MICs = 30 and 15  $\mu$ M, respectively).<sup>2</sup> It is cytotoxic to HepG2 cells (IC<sub>50</sub> = 18  $\mu$ M).

### References

1. Bols, M., Biderup, L., Hansen, J., *et al.* Inhibition of collagenase by aranciamycin and aranciamycin derivatives. *J. Med. Chem.* **35**(15), 1768-2771 (1992).
2. 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), 949-52 (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, 12/13/2022

#### 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