

# PRODUCT INFORMATION



## Anguinomycin A

Item No. 20598

CAS Registry No.: 111278-01-4  
Formal Name: 19-(3,6-dihydro-6-oxo-2H-pyran-2-yl)-6-hydroxy-3,5,7,9,11,15,17-heptamethyl-8-oxo-2,10,12,16,18-nonadecapentaenoic acid

Synonym: 5-Demethylleptomycin A

MF: C<sub>31</sub>H<sub>44</sub>O<sub>6</sub>

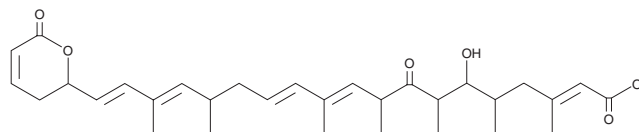
FW: 512.7

Purity: ≥95%

Supplied as: A film

Storage: -20°C

Stability: ≥4 years



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

### Laboratory Procedures

Anguinomycin A is supplied as a film. A stock solution may be made by dissolving the anguinomycin A in the solvent of choice, which should be purged with an inert gas. Anguinomycin A is soluble in organic solvents such as ethanol and methanol.

### Description

Anguinomycin A is an antibiotic first isolated from a *Streptomyces* sp. and analog of leptomycin B (Item No. 10004976) that is highly cytotoxic to mouse P388 leukemia cells *in vitro* (IC<sub>50</sub> = 0.1-0.2 ng/ml).<sup>1</sup> It also demonstrates antitumor activity at 62.5 µg/kg/day against Lewis lung carcinoma in mice.<sup>1</sup>

### References

1. Hayakawa, Y., Adachi, K., and Komeshima, N. New antitumor antibiotics, anguinomycins A and B. *J. Antibiot. (Tokyo)* **40(9)**, 1349-1352 (1987).

#### 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, 11/11/2022

#### 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