

# PRODUCT INFORMATION

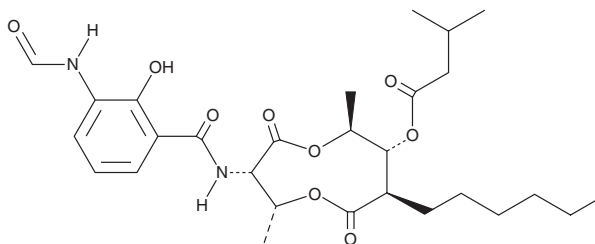


## Antimycin A Complex

Item No. 34799

**CAS Registry No.:** 1397-94-0  
**Formal Name:** (2R,3S,6S,7R,8R)-3-(3-formamido-2-hydroxybenzamido)-8-hexyl-2,6-dimethyl-4,9-dioxo-1,5-dioxonan-7-yl 3-methylbutanoate

**MF:** C<sub>28</sub>H<sub>40</sub>N<sub>2</sub>O<sub>9</sub> (for Antimycin A<sub>1</sub>)  
**FW:** 548.6  
**Purity:** ≥95% (mixture of subunits)  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥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

Antimycin A complex is supplied as a solid. A stock solution may be made by dissolving the antimycin A complex in the solvent of choice, which should be purged with an inert gas. Antimycin A complex is soluble in the organic solvent ethanol.

### Description

Antimycin A complex is a bacterial metabolite complex that has been found in *Streptomyces* and is composed of the major components antimycin A<sub>1</sub> (Item No. 19433), -A<sub>2</sub> (Item No. 21997), -A<sub>3</sub> (Item No. 21998), and -A<sub>4</sub> (Item No. 21999), as well as minor components, including antimycin A<sub>5</sub> and -A<sub>6</sub>, among others.<sup>1</sup> It induces the synthesis of carotenoids in *M. marinum* when used at concentrations ranging from 3.3 to 160 μM.<sup>2</sup> Antimycin A complex (25-125 ppb) has teleocidal activity against goldfish.<sup>1</sup>

### References

1. Kluepfel, D., Sehgal, S.N., and Vezina, C. Antimycin A components. I. Isolation and biological activity. *J. Antibiot. (Tokyo)* **23(2)**, 75-80 (1970).
2. Batra, P.P., Harbin, T.L., Howes, C.D., et al. A study of the relationship of structure and activity of antimycin A in the induction of carotenoid synthesis in *Mycobacterium marinum*. *J. Biol. Chem.* **246(23)**, 7125-7130 (1971).

#### 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, 12/15/2022

#### 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