

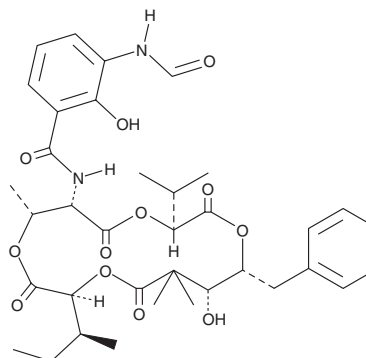
# PRODUCT INFORMATION



## Neoantimycin

Item No. 23645

**CAS Registry No.:** 22862-63-1  
**Formal Name:** 3-(formylamino)-2-hydroxy-N-[(3S,6S,7R,10S,14R,15R)-14-hydroxy-7,13,13-trimethyl-3-(1-methylethyl)-10-[(1S)-1-methylpropyl]-2,5,9,12-tetraoxo-15-(phenylmethyl)-1,4,8,11-tetraoxacyclopentadec-6-yl]-benzamide  
**MF:**  $C_{36}H_{46}N_2O_{12}$   
**FW:** 698.8  
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**Storage:**  $-20^{\circ}\text{C}$   
**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

Neoantimycin is supplied as a solid. A stock solution may be made by dissolving the neoantimycin in the solvent of choice, which should be purged with an inert gas. Neoantimycin is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Neoantimycin is an antibiotic originally isolated from *S. orinoci* that has slight antifungal activity.<sup>1</sup> It is structurally similar to antimycin (Item No. 19433) despite being produced by an unrelated species of *Streptomyces*. The biosynthetic pathway of neoantimycin has been determined and active analogs discovered.<sup>2,3</sup>

### References

1. Cassinelli, G., Grein, A., Orezzi, P., et al. New antibiotics produced by *Streptoverticillium orinoci*, n. sp. *Arch. Mikrobiol.* **55**(4), 358-368 (1967).
2. Li, X., Zvanych, R., Vanner, S.A., et al. Chemical variation from the neoantimycin depsipeptide assembly line. *Bioorg. Med. Chem. Lett* **23**(18), 5123-5127 (2013).
3. Umeda, Y., Chijiwa, S., Furihata, K., et al. Prunustatin A, a novel GRP78 molecular chaperone down-regulator isolated from *Streptomyces violaceoniger*. *J. Antibiot. (Tokyo)* **58**(3), 206-209 (2005).

#### 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, 01/03/2023

#### 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](http://WWW.CAYMANCHEM.COM)