

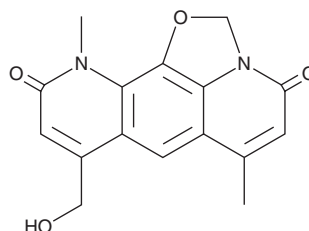
# PRODUCT INFORMATION



## Nybomycin

Item No. 28099

CAS Registry No.: 30408-30-1  
Formal Name: 8-(hydroxymethyl)-6,11-dimethyl-2H,4H-oxazolo[5,4,3-ij]pyrido[3,2-g]quinoline-4,10(11H)-dione  
Synonym: NSC 613948  
MF:  $C_{16}H_{14}N_2O_4$   
FW: 298.3  
Purity:  $\geq 95\%$   
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

Nybomycin is supplied as a solid. A stock solution may be made by dissolving the nybomycin in the solvent of choice, which should be purged with an inert gas. Nybomycin is soluble in the organic solvent acetic acid (warmed).

### Description

Nybomycin is a fungal metabolite originally isolated from *Streptomyces* sp. A717 and has antibacterial activity.<sup>1</sup> It is active against *B. subtilis*, *B. cereus*, *B. mycoides*, *M. smegmatis*, and *K. pneumoniae* (MICs = 0.025, 2, 0.8, 0.2, and 0.003  $\mu\text{g/ml}$ , respectively). Nybomycin is also active against *M. tuberculosis* and *M. bovis* (MICs = 4.5 and 1  $\mu\text{g/ml}$ , respectively).<sup>2</sup>

### References

1. Strelitz, F., Flon, H., and Asheshov, I.N. Nybomycin, a new antibiotic with antiphage and antibacterial properties. *Proc. Natl. Acad. Sci. U.S.A.* **41**(9), 620-624 (1955).
2. Arai, M., Kamiya, K., Pruksakorn, P., et al. Anti-dormant mycobacterial activity and target analysis of nybomycin produced by a marine-derived *Streptomyces* sp. *Bioorg. Med. Chem.* **23**(13), 3534-3541 (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, 11/03/2022

#### 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