

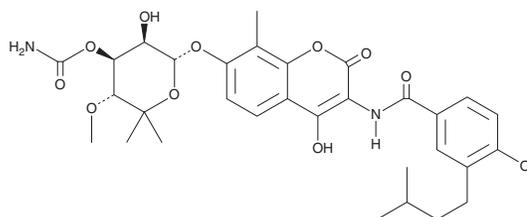
PRODUCT INFORMATION



Dihydronovobiocin

Item No. 29649

CAS Registry No.: 29826-16-2
Formal Name: N-[7-[[3-O-(aminocarbonyl)-6-deoxy-5-C-methyl-4-O-methyl- α -L-lyxo-hexopyranosyl]oxy]-4-hydroxy-8-methyl-2-oxo-2H-1-benzopyran-3-yl]-4-hydroxy-3-(3-methylbutyl)-benzamide



MF: C₃₁H₃₈N₂O₁₁
FW: 614.6
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years

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

Laboratory Procedures

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

Description

Dihydronovobiocin is a coumarin antibiotic and derivative of novobiocin (Item No. 18457).¹ It is active against the bacteria *S. aureus*, *S. haemolyticus*, *D. pneumoniae*, *S. typhosa*, *K. pneumoniae*, and *P. multocida* (MICs = 0.6, 2, 0.6, 10, 10, and 3 μ g/ml, respectively).² Dihydronovobiocin inhibits DNA gyrase subunit B with an IC₅₀ value of 64.5 nM.³

References

- Berger, J. and Batcho, A.D. Coumarin - glycoside antibiotics. *J. Chromatogr. Lib.* **15**, 101-158 (1978).
- Hoeksema, H. Dihydronovobiocin and derivatives thereof. *The Upjohn Company. US3175944*, (1965).
- Gevi, M. and Domenici, E. A scintillation proximity assay amenable for screening and characterization of DNA gyrase B subunit inhibitors. *Anal. Biochem.* **300(1)**, 34-39 (2002).

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, 10/13/2022

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