

PRODUCT INFORMATION



Nodusmicin

Item No. 23147

CAS Registry No.: 76265-48-0
Formal Name: (1E,3R,4S,7S,8aS,10aR,11R,12R,13R,14R,14aS,14bS)-3,4,8,8a,10a,11,12,13,14,14a-decahydro-12,14-dihydroxy-4-[(1R)-1-hydroxyethyl]-7-methoxy-1,3,13-trimethyl-11,14b-epoxy-14bH-naphth[2,1-e]oxecin-6(7H)-one

MF: C₂₃H₃₄O₇

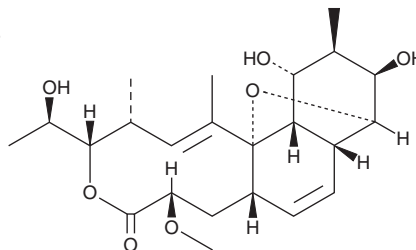
FW: 422.5

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

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

Description

Nodusmicin is a macrolide antibiotic.¹ It has activity against antibiotic-susceptible and -resistant *S. aureus* with MIC values of 125, 250, and 250 µg/ml for UC-76, UC-6685, and UC-6690 strains, respectively.

Reference

1. Magerlein, B.J. and Mizsak, S.A. 9-O-Esters of nodusmicin. *J. Antibiot.* **35**(1), 111-112 (1982).

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, 09/30/2022

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