

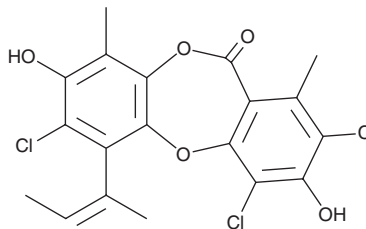
# PRODUCT INFORMATION



## Nornidulin

Item No. 23646

CAS Registry No.: 33403-37-1  
Formal Name: 2,4,7-trichloro-3,8-dihydroxy-1,9-dimethyl-6-(1-methyl-1-propen-1-yl)-11H-dibenzo[b,e][1,4]dioxepin-11-one  
MF:  $C_{19}H_{15}Cl_3O_5$   
FW: 429.7  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years  
Item Origin: Fungus/*Emericella* sp.



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

### Laboratory Procedures

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

### Description

Nornidulin is a depsidone originally isolated from *A. nidulans* that has antibacterial activity against *M. tuberculosis* and *M. ranee* as well as antifungal activity against *T. tonsurans* and *M. audouini*.<sup>1</sup> It also inhibits the growth of methicillin-resistant *S. aureus* (MRSA; MIC = 2  $\mu g/ml$ ).<sup>2</sup> Nornidulin has cytotoxic activity in MOLT-3 cells (IC<sub>50</sub> = 35.7  $\mu M$ ) but not HuCCA-1, HepG2, or A549 cells (IC<sub>50</sub>s = >116.4  $\mu M$ ).<sup>3</sup>

### References

1. Dean, F.M., Roberston, A., Roberts, J.C., *et al.* Nidulin and 'Ustin': Two chlorine-containing metabolic products of *Aspergillus nidulans*. *Nature* **172**(4372), 344 (1953).
2. Zhang, Y., Mu, J., Feng, Y., *et al.* Four chlorinated depsidones from a seaweed-derived strain of *Aspergillus unguis* and their new biological activities. *Nat. Prod. Res.* **28**(7), 503-506 (2014).
3. Sureram, S., Wiyakrutta, S., Ngamrojanavanich, N., *et al.* Depsidones, aromatase inhibitors and radical scavenging agents from the marine-derived fungus *Aspergillus unguis* CRI282-03. *Planta Med.* **78**(6), 582-588 (2012).

#### 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, 08/25/2023

#### 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