

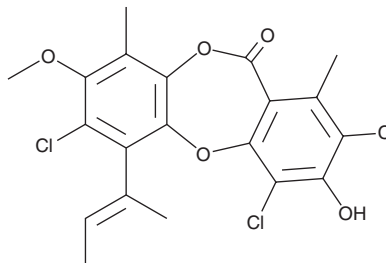
PRODUCT INFORMATION



Nidulin

Item No. 27955

CAS Registry No.: 10089-10-8
Formal Name: 2,4,7-trichloro-3-hydroxy-8-methoxy-1,9-dimethyl-6-[(1E)-1-methyl-1-propen-1-yl]-11H-dibenzo[b,e][1,4]dioxepin-11-one
Synonym: Methylustin
MF: C₂₀H₁₇Cl₃O₅
FW: 443.7
Purity: ≥99%
Supplied as: A solid
Storage: -20°C
Stability: ≥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

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

Nidulin is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Nidulin is a depsidone originally isolated from *A. nidulans*.¹ It is active against the bacteria *M. tuberculosis* and *M. ranoe*, as well as the fungi *T. tonsurans* and *M. audouini*. It also inhibits the growth of methicillin-resistant *S. aureus* (MRSA; MIC = 4 µg/ml).² Nidulin is cytotoxic to MOLT-3 cells (IC₅₀ = 21.2 µM) but not HuCCA-1, HepG2, or A549 cells (IC₅₀s = >112.7 µM). It inhibits aromatase with an IC₅₀ value of 11.2 µM.³

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, 08/28/2023

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