

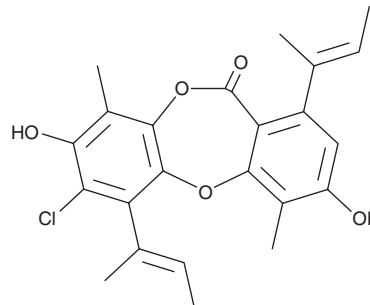
PRODUCT INFORMATION



Emeguisin A

Item No. 32929

CAS Registry No.: 117032-54-9
Formal Name: 7-chloro-3,8-dihydroxy-4,9-dimethyl-1,6-bis(1-methyl-1-propenyl)-11H-dibenzo[b,e][1,4]dioxepin-11-one
MF: C₂₃H₂₃ClO₅
FW: 414.9
Purity: ≥95%
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

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

Description

Emeguisin A is a depsidone fungal metabolite originally isolated from *E. unguis*.¹ It is active against the bacteria *S. aureus* and methicillin-resistant *S. aureus* (MRSA; MIC = 0.5 µg/ml for both), the fungus *C. neoformans* (MIC = 0.5 µg/ml), and the protozoan *P. falciparum* (IC₅₀ = 2.2 µM).²

References

1. Kawahara, N., Nozawa, K., Nakajima, S., *et al.* Isolation and structures of novel fungal depsidones, emeguisins A, B, and C, from *Emericella unguis*. *J. Chem. Soc., Perkin Trans. 1* **9**, 2611-2614 (1988).
2. Klaiklay, S., Rukachaisirikul, V., Aungphao, W., *et al.* Depsidone and phthalide derivatives from the soil-derived fungus *Aspergillus unguis* PSU-RSPG199. *Tetrahedron Lett.* **57(39)**, 4348-4351 (2016).

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, 08/17/2022

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