# PRODUCT INFORMATION



## **Unguinol**

Item No. 24207

CAS Registry No.: 36587-59-4

Formal Name: 3,8-dihydroxy-1,9-dimethyl-6-[(1E)-1-

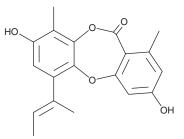
methyl-1-propen-1-yl]-11H-dibenzo[b,e]

[1,4]dioxepin-11-one

MF:  $C_{19}H_{18}O_{5}$ FW: 326.3 **Purity:** ≥95% Supplied as: A solid Storage: -20°C Stability: ≥4 years

Item Origin: Fungus/Aspergillus sp.

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



#### **Laboratory Procedures**

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

#### Description

Unguinol is a depsidone originally isolated from A. unguis. 1 It is an inhibitor of pyruvate phosphate dikinase (PPDK;  $IC_{50} = 42.3 \mu M$ ).<sup>2</sup> It inhibits the growth of plants utilizing  $C_4$ , but not  $C_3$ , carbon fixation. Unguinol also inhibits the growth of the bacteria *S. aureus* and *V. harveyi* ( $GI_{50}s = 8.7$  and  $69.5 \mu M$ , respectively) and H460, MCF-7, and SF-268 cancer cells ( $GI_{50}s = 28.2$ , 50.8, and 44.3  $\mu M$ , respectively).

#### References

- 1. Stodola, F.H., Vesonder, R.F., Fennell, D.I., et al. New depsidone from Aspergillus unguis. Phytochemistry **11(6)**, 2107-2108 (1972).
- 2. Motti, C.A., Bourne, D.G., Burnell, J.N., et al. Screening marine fungi for inhibitors of the C<sub>4</sub> plant enzyme pyruvate phosphate dikinase: Unguinol as a potential novel herbicide candidate. Appl. Environ. Microbiol. 73(6), 1921-1927 (2007).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information Buyer agrees to purchase the material can be found on our website.

Copyright Cayman Chemical Company, 08/10/2023

### **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