

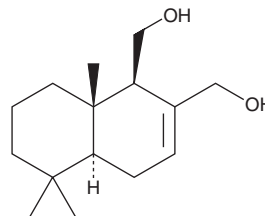
PRODUCT INFORMATION



Drimendiol

Item No. 29820

CAS Registry No.: 34437-62-2
Formal Name: (1R,4aS,8aS)-1,4,4a,5,6,7,8,8a-octahydro-5,5,8a-trimethyl-1,2-naphthalenedimethanol
MF: C₁₅H₂₆O₂
FW: 238.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Semi-synthetic



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

Laboratory Procedures

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

Description

Drimendiol is a sesquiterpene that has been found in *W. ugandensis*.¹ It inhibits *C. albicans*, *S. aureus*, and *S. epidermidis* biofilm formation with 50% biofilm inhibitory concentration values (BIC_{50S}) of 25.5, 65.1, and 67.1 µg/ml, respectively. It has antifouling activity, inhibiting the settlement of *C. savignyi* and *B. improvisus* larvae on a Petri dish surface (EC_{50S} = 1 and 0.5 µg/ml, respectively).²

References

1. Kipanga, P.N., Liu, M., Panda, S.K., et al. Biofilm inhibiting properties of compounds from the leaves of *Warburgia ugandensis* Sprague subsp *ugandensis* against *Candida* and staphylococcal biofilms. *J. Ethnopharmacol.* **248:112352** (2020).
2. Moodie, L.W., Trepos, R., Cervin, G., et al. Probing the structure-activity relationship of the natural antifouling agent polygodial against both micro- and macrofoulers by semisynthetic modification. *J. Nat. Prod.* **80(2)**, 515-525 (2017).

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, 10/21/2022

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