

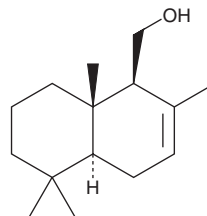
# PRODUCT INFORMATION



## (-)-Drimenol

Item No. 29428

**CAS Registry No.:** 468-68-8  
**Formal Name:** (1S,4aS,8aS)-1,4,4a,5,6,7,8,8a-octahydro-2,5,5,8a-tetramethyl-1-naphthalenemethanol  
**Synonym:** NSC 169775  
**MF:** C<sub>15</sub>H<sub>26</sub>O  
**FW:** 222.4  
**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

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

### Description

(-)-Drimenol is a sesquiterpene alcohol originally isolated from *B. trilobata* that has antimicrobial and anticancer activities.<sup>1-3</sup> It is active against the bacteria *S. aureus*, *B. cereus*, *A. baumannii*, *E. coli*, and *P. aeruginosa* (MICs = 667, 667, 583, 1,333, and 667 µg/ml, respectively), as well as the fungi *C. cucumerinum* and *S. tritici* (IC<sub>50</sub>s = 6.6 and 80.1 µg/ml, respectively).<sup>1,2</sup> (-)-Drimenol decreases the proliferation of A2058 and A375 melanoma cell lines with IC<sub>50</sub> values of 33.5 and 31.25 µg/ml, respectively.<sup>3</sup>

### References

1. Santos, T.G., Dognnini, J., Begnini, I.M., *et al.* Chemical characterization of essential oils from *Drimys angustifolia* miers (winteraceae) and antibacterial activity of their major compounds. *J. Braz. Chem. Soc.* **24(1)**, 164-170 (2013).
2. Scher, J.M., Speakman, J.B., Zapp, J., *et al.* Bioactivity guided isolation of antifungal compounds from the liverwort *Bazzania trilobata* (L.) S.F. Gray. *Phytochemistry* **65(18)**, 2583-2588 (2004).
3. Russo, A., Cardile, V., Graziano, A.C.E., *et al.* Antigrowth activity and induction of apoptosis in human melanoma cells by *Drymis winteri* forst extract and its active components. *Chem. Biol. Interact.* **305**, 79-85 (2019).

#### 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, 12/06/2022

#### 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