

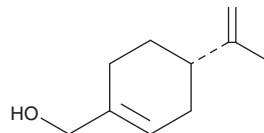
# PRODUCT INFORMATION



## (-)-Perillyl Alcohol

Item No. 21860

<b>CAS Registry No.:</b>	18457-55-1
<b>Formal Name:</b>	4S-(1-methylethenyl)-1-cyclohexene-1-methanol
<b>Synonyms:</b>	(L)-Perillyl Alcohol, (S)-Perillic Alcohol, (S)-(-)-Perillyl Alcohol
<b>MF:</b>	C <sub>10</sub> H <sub>16</sub> O
<b>FW:</b>	152.2
<b>Purity:</b>	≥95%
<b>Supplied as:</b>	A neat oil
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥4 years
<b>Item Origin:</b>	Synthetic



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

### Laboratory Procedures

(-)-Perillyl alcohol is supplied as a neat oil. A stock solution may be made by dissolving the (-)-perillyl alcohol in the solvent of choice, which should be purged with an inert gas. (-)-Perillyl alcohol is miscible in organic solvents such as ethanol, DMSO, and dimethyl formamide.

### Description

(-)-Perillyl alcohol is a monoterpene alcohol that has been found in lavender essential oil and has diverse biological activities.<sup>1-4</sup> It reduces production of hydroperoxidiene and thiobarbituric acid reactive substances (TBARS) *in vitro* in a concentration-dependent manner.<sup>2</sup> (-)-Perillyl alcohol inhibits PANC-1 pancreatic carcinoma cell and H-Ras-transformed fibroblast growth when used at a concentration of 1 mM.<sup>3</sup> It also inhibits the growth of *P. aeruginosa*, *E. coli*, *S. aureus*, and *C. albicans* (MICs = 480-2,900 ppm).<sup>4</sup>

### References

- Gerhäuser, C., Klimo, K., Heiss, E., *et al.* Mechanism-based *in vitro* screening of potential cancer chemopreventive agents. *Mutat. Res.* **523-524**, 163-172 (2003).
- Ruberto, G. and Baratta, M.T. Antioxidant activity of selected essential oil components in two lipid model systems. *Food Chem.* **69(2)**, 167-174 (2000).
- Karlson, J., Borg-Karlson, A.K., Unelius, R., *et al.* Inhibition of tumor cell growth by monoterpenes *in vitro*: Evidence of a Ras-independent mechanism of action. *Anticancer Drugs* **7(4)**, 422-429 (1996).
- Griffin, S.G., Wyllie, S.G., Markham, J.L., *et al.* The role of structure and molecular properties of terpenoids in determining their antimicrobial activity. *Flavour Fragr. J.* **14(5)**, 322-332 (1999).

#### 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, 11/14/2022

#### 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