

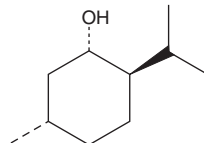
PRODUCT INFORMATION



(+)-Menthol

Item No. 23463

CAS Registry No.: 15356-60-2
Formal Name: 5S-methyl-2R-(1S-methylethyl)-cyclohexanol
Synonym: D-Menthol
MF: C₁₀H₂₀O
FW: 156.3
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

(+)-Menthol is supplied as a crystalline solid. A stock solution may be made by dissolving the (+)-menthol in the solvent of choice. (+)-Menthol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of (+)-menthol in these solvents is approximately 30 mg/ml.

Description

(+)-Menthol is a monoterpene alcohol that has been found in *Cannabis* and has antifungal activity.^{1,2} It inhibits the growth of *F. verticillioides* (MIC = 1.5 mM).² Unlike (-)-menthol, (+)-menthol does not exhibit analgesia, antibacterial, anticancer, or cholinesterase inhibitory activities.³

References

1. Elzinga, S., Fishedick, J., Podkolinski, R., *et al.* Cannabinoids and terpenes as chemotaxonomic markers in cannabis. *Nat. Prod. Chem. Res.* **3(4)**, 181 (2015).
2. Dambolena, J.S., López, A.G., Rubinstein, H.R., *et al.* Effects of menthol stereoisomers on the growth, sporulation and fumonisin B1 production of *Fusarium verticillioides*. *Food Chem.* **123(1)**, 165-170 (2010).
3. Kamatou, G.P., Vermaak, I., Viljoen, A.M., *et al.* Menthol: A simple monoterpene with remarkable biological properties. *Phytochemistry* **96**, 15-25 (2013).

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, 10/03/2022

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