

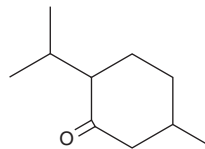
PRODUCT INFORMATION



Menthone

Item No. 34683

CAS Registry No.: 10458-14-7
Formal Name: 5-methyl-2-(1-methylethyl)-cyclohexanone
Synonyms: *p*-Menthan-3-one, NSC 113134, NSC 9280
MF: C₁₀H₁₈O
FW: 154.3
Purity: ≥95%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years
Item Origin: Synthetic



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

Laboratory Procedures

Menthone is supplied as a liquid. A stock solution may be made by dissolving the menthone in the solvent of choice, which should be purged with an inert gas. Menthone is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

Menthone is a monoterpene that has been found in *A. rugosa* and has diverse biological activities.¹⁻⁴ It inhibits the seed germination of *R. sativus* when used at concentrations of 0.1 and 1 mM.² Menthone has insecticidal activity against *C. molestus* fourth-instar larvae (LC₅₀ = 156 mg/L).³ It stimulates the proliferation of isolated human T lymphocytes derived from allergic patients when used at concentrations of 6.2 and 62 µg/ml.⁴

References

1. Haiyan, G., Lijuan, H., Shaoyu, L., *et al.* Antimicrobial, antibiofilm and antitumor activities of essential oil of *Agastache rugosa* from Xinjiang, China. *Saudi J. Biol. Sci.* **23(4)**, 524-530 (2016).
2. De Martino, L., Mancini, E., de Almeida, L.F., *et al.* The antigerminative activity of twenty-seven monoterpenes. *Molecules* **15(9)**, 6630-6637 (2010).
3. Traboulsi, A.F., Taoubi, K., el-Haj, S., *et al.* Insecticidal properties of essential plant oils against the mosquito *Culex pipiens molestus* (Diptera: Culicidae). *Pest Manag. Sci.* **58(5)**, 491-495 (2002).
4. Cariddi, L., Escobar, F., Moser, M., *et al.* Monoterpenes isolated from *Minthostachys verticillata* (Griseb.) Epling essential oil modulates immediate-type hypersensitivity responses *in vitro* and *in vivo*. *Planta Med.* **77(15)**, 1687-1694 (2011).

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, 08/24/2021

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