

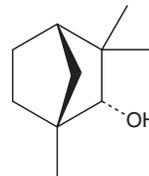
PRODUCT INFORMATION



(+)-Fenchol

Item No. 23162

CAS Registry No.: 2217-02-9
Formal Name: (1R,2R,4S)-1,3,3-trimethyl-bicyclo[2.2.1]heptan-2-ol
Synonyms: endo-(+)-Fenchyl Alcohol, (1R)-endo-(+)-Fenchyl Alcohol, (+)- α -Fenchol, D-Fenchyl Alcohol, 1,3,3-Trimethyl-2-norbornanol
MF: C₁₀H₁₈O
FW: 154.3
Purity: \geq 95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

(+)-Fenchol is supplied as a crystalline solid. A stock solution may be made by dissolving the (+)-fenchol in the solvent of choice, which should be purged with an inert gas. (+)-Fenchol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (+)-fenchol in ethanol and DMSO is approximately 16 mg/ml and approximately 25 mg/ml in DMF.

(+)-Fenchol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (+)-fenchol should first be dissolved in DMF and then diluted with the aqueous buffer of choice. (+)-Fenchol has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(+)-Fenchol is a monoterpene that has been found in plants, including *Cannabis*.^{1,2} It has been used in the synthesis of other terpenoids and as a chiral building block in organic synthesis of various compounds.^{3,4} Formulations containing (+)-fenchol have been used as fragrance ingredients.

Reference

1. Mothana, R.A., Al-Said, M.S., Al-Yahya, M.A., *et al.* GC and GC/MS analysis of essential oil composition of the endemic Soqatraen *Leucas virgata* Balf.f. and its antimicrobial and antioxidant activities. *Int. J. Mol. Sci.* **14**(11), 23129-23139 (2013).
2. Elzinga, S., Fishedick, J., Podkolinski, R., *et al.* Cannabinoids and terpenes as chemotaxonomic markers in cannabis. *Nat. Prod. Chem. Res.* **3**(4), 181 (2015).
3. van der Zeijden, A.A.H. and Mattheis, C. Synthesis of chiral O-functionalized isobornyloxy, menthyloxy and fenchyloxy cyclopentadienyl ligands. *Synthesis* **1996**(7), 847-850 (1996).
4. Yuasa, Y., Watanabe, T., Nagakura, A., *et al.* An improved synthesis of (S)-aspartyl-(7,7-dimethylnorborn-2R-yl)-(S)-alanine methyl ester, a new high intensity artificial sweetener. *Tetrahedron* **48**(17), 3473-3484 (1992).

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, 11/11/2022

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