PRODUCT INFORMATION



(+)-β-Citronellol

Item No. 23464

CAS Registry No.:	1117-61-9
Formal Name:	3R,7-dimethyl-6-octen-1-ol
Synonyms:	(R)-Citronellol, (+)-Citronellol,
	(+)-(R)-Citronellol, (R)-(+)-β-Citronellol
MF:	C ₁₀ H ₂₀ O
FW:	156.3
Purity:	≥98%
Supplied as:	A liquid
Storage:	-20°C
Stability:	≥2 years
Information represent	the product specifications. Batch specific analytical results are provided on each certificate of analysi

Laboratory Procedures

(+)- β -Citronellol is supplied as a liquid. A stock solution may be made by dissolving the (+)- β -citronellol in the solvent of choice, which should be purged with an inert gas. (+)- β -Citronellol is soluble in chloroform.

(+)-β-Citronellol is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

(+)-β-Citronellol is a monoterpene that has been found in Cannabis.¹ It inhibits degranulation of cultured mast cells by 21.3% when used at a concentration of 0.5 mM.² Unlike (R)-citronellal, it does not reduce the perceived bitterness of caffeine when used at a concentration of 5.9 nM.³ Formulations containing (+)- β -citronellol have been used as fragrance ingredients in various cosmetics, toiletries, and cleaning products.

References

- 1. Fletcher, R.S. and McKay, J. Industrial hemp cannabis cultivars and seeds with stable cannabinoid profiles. New West Genetics US2017/0339907A1 (2017).
- 2. Kobayashi, Y., Sato, H., Yorita, M., et al. Inhibitory effects of geranium essential oil and its major component, citronellol, on degranulation and cytokine production by mast cells. Biosci. Biotechnol. Biochem. 80(6), 1172-1178 (2016).
- 3. Suess, B., Brockhoff, A., Meyerhof, W., et al. The odorant (R)-citronellal attenuates caffeine bitterness by inhibiting the bitter receptors TAS2R43 and TAS2R46. J. Agric. Food Chem. 66(10), 2301-2311 (2018).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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

uyer 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, 11/08/2022

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