PRODUCT INFORMATION



Nerol

Item No. 25774

CAS Registry No.:	106-25-2
Formal Name:	3,7-dimethyl-2Z,6-octadien-1-ol
Synonyms:	(Z)-Geraniol, <i>cis</i> -Geraniol, β-Nerol
MF:	C ₁₀ H ₁₈ O
FW:	154.3
Purity:	≥95%
Supplied as:	An oil
Storage:	4°C
Stability:	≥4 years
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.	

Laboratory Procedures

Nerol is supplied as an oil. A stock solution may be made by dissolving the nerol in the solvent of choice. Nerol is slightly soluble in chloroform and methanol.

Description

Nerol is a monoterpene and isomer of geraniol (Item No. 23166) that has been found in a variety of plants, including Cannabis.¹ It increases production of reactive oxygen species (ROS) and intracellular calcium levels and induces mitochondrial dysfunction, cytochrome C release, and apoptosis in A. flavus.² Nerol (30-300 mg/kg) reduces weight loss, production of the inflammatory cytokines IL-13 and TNF- α , gastric damage, and hyperalgesia in a mouse model of oxazolone-induced colitis.³

References

- 1. Elzinga, S., Fischedick, J., Podkolinski, R., et al. Cannabinoids and terpenes as chemotaxonomic markers in Cannabis. Nat. Prod. Chem. Res. 3(4), 181 (2015).
- 2. Tian, J., Gan, Y., Pan, C., et al. Nerol-induced apoptosis associated with the generation of ROS and Ca²⁺ overload in saprotrophic fungus Aspergillus flavus. Appl. Microbiol. Biotechnol. 102(15), 6659-6672 (2018).
- 3. González-Ramírez, A.E., González-Trujano, M.E., Orozco-Suárez, S.A., et al. Nerol alleviates pathologic markers in the oxazolone-induced colitis model. Eur. J. Pharmacol. 776, 81-89 (2016).

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.

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, 11/08/2022

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