

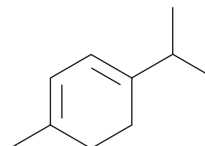
PRODUCT INFORMATION



α -Terpinene

Item No. 23178

CAS Registry No.: 99-86-5
Formal Name: 1-methyl-4-(1-methylethyl)-1,3-cyclohexadiene
Synonym: *p*-Mentha-1,3-diene
MF: C₁₀H₁₆
FW: 136.2
Purity: $\geq 95\%$
Supplied as: A liquid
Storage: -20°C
Stability: ≥ 2 years



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

Laboratory Procedures

α -Terpinene is supplied as a liquid. A stock solution may be made by dissolving the α -terpinene in the solvent of choice, which should be purged with an inert gas. α -Terpinene is soluble in the organic solvent ethanol at a concentration of approximately 100 mg/ml.

Description

α -Terpinene is a terpenoid that has been found in *Cannabis* and has diverse biological activities, including acaricidal, antiprotozoal, and antioxidant properties.^{1,2} α -Terpinene induces 99 and 46% mortality of male and female *T. putrescentiae* larvae, respectively, when administered at a dose of 32.3 μ L/L via inhalation.³ In mice infected with *T. evansi*, it increases longevity when used alone and decreases mortality when used in combination with diminazene aceturate (Item No. 18678).⁴ It also scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals when used at a concentration of 10 mM.⁵

References

1. Giese, M.W., Lewis, M.A., Giese, L., *et al.* Development and validation of a reliable and robust method for the analysis of cannabinoids and terpenes in Cannabis. *J. AOAC Int.* **98(6)**, 1503-1522 (2015).
2. Ross, S.A. and ElSohly, M.A. The volatile oil composition of fresh and air-dried buds of *Cannabis sativa*. *J. Nat. Prod.* **59(1)**, 49-51 (1996).
3. Sánchez-Ramos, I. and Castañera, P. Acaricidal activity of natural monoterpenes on *Tyrophagus putrescentiae* (Schrank), a mite of stored food. *J. Stored Prod. Res.* **37(1)**, 93-101 (2000).
4. Baldissera, M.D., Grando, T.H., Souza, C.F., *et al.* *In vitro* and *in vivo* action of terpinen-4-ol, γ -terpinene, and α -terpinene against *Trypanosoma evansi*. *Exp. Parasitol.* **162**, 43-48 (2016).
5. Kim, H.J., Chen, F., Wu, C., *et al.* Evaluation of antioxidant activity of Australian tea tree (*Melaleuca alternifolia*) oil and its components. *J. Agric. Food Chem.* **52(10)**, 2849-2854 (2004).

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, 11/09/2022

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