

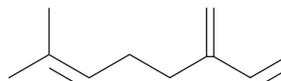
PRODUCT INFORMATION



β-Myrcene

Item No. 21720

CAS Registry No.: 123-35-3
Formal Name: 7-methyl-3-methylene-1,6-octadiene
Synonym: NSC 406264
MF: C₁₀H₁₆
FW: 136.2
Purity: ≥95%
Supplied as: A liquid
Storage: 4°C
Stability: ≥2 years



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

Laboratory Procedures

β-Myrcene is supplied as a liquid. A stock solution may be made by diluting the β-myrcene in the solvent of choice, which should be purged with an inert gas. β-Myrcene is miscible in organic solvents such as ethanol, DMSO, and chloroform.

Description

β-Myrcene is a terpene that has been found in *Cannabis* and has antioxidative properties.¹ It reduces thiobarbituric acid reactive substance (TBARS) and increases glutathione (GSH), catalase, GSH peroxidase, and CuZn superoxide dismutase levels in a rat model of persistent environmental pollutant-induced oxidative stress when administered at a dose of 200 mg/kg per day.²

References

1. Lewis, M.A., Russo, E.B., and Smith, K.M. Pharmacological foundations of Cannabis chemovars. *Planta Med.* **84**(4), 225-233 (2018).
2. Ciftci, O., Ozdemir, I., Tanyildizi, S., *et al.* Antioxidative effects of curcumin, β-myrcene and 1,8-cineole against 2,3,7,8-tetrachlorodibenzo-*p*-dioxin-induced oxidative stress in rats liver. *Toxicol. Ind. Health* **27**(5), 447-453 (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, 03/12/2020

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