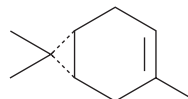


PRODUCT INFORMATION

(+)-3-Carene

Item No. 23165

CAS Registry No.: 498-15-7
 Formal Name: (1S,6R)-3,7,7-trimethyl-bicyclo[4.1.0]hept-3-ene
 Synonym: (S)-(+)-3-Carene
 MF: C₁₀H₁₆
 FW: 136.2
 Purity: ≥98%
 Supplied as: A neat oil
 Storage: -20°C
 Stability: ≥4 years



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

Laboratory Procedures

(+)-3-Carene is supplied as a neat oil. A stock solution may be made by dissolving the (+)-3-carene in the solvent of choice, which should be purged with an inert gas. (+)-3-Carene is soluble in organic solvents such as chloroform and DMSO. The solubility of (+)-3-carene in these solvents is approximately 10 mg/ml.

Description

(+)-3-Carene is a monoterpene that has been found in various plants, including *C. sativa*, *C. indica*, and hemp.¹ It is a component of (±)-3-carene (Item No. 25768).

Reference

1. Hazekamp, A., Tejkalová, K., and Papadimitriou, S. *Cannabis*: From cultivar to chemovar II—A metabolomics approach to *Cannabis* classification. *Cannabis Cannabinoid Res.* **1**(1), 202-215 (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.

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, 09/14/2023

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