

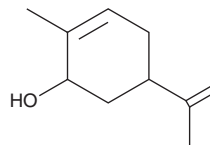
PRODUCT INFORMATION



Carveol

Item No. 33831

CAS Registry No.: 99-48-9
Formal Name: 2-methyl-5-(1-methylethenyl)-2-cyclohexen-1-ol
Synonym: NSC 68313
MF: C₁₀H₁₆O
FW: 152.2
Purity: ≥98%
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

Carveol is supplied as a liquid. A stock solution may be made by dissolving the carveol in the solvent of choice, which should be purged with an inert gas. Carveol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of carveol in these solvents is approximately 30 mg/ml.

Description

Carveol is a monoterpene alcohol with nematocidal activity.¹ It reduces hatching in *M. incognita* egg masses and reduces mobility of *M. incognita* second-stage juveniles when used at concentrations of 250 and 500 mg/L. Soil application of carveol (250 mg/kg) reduces *M. incognita*-induced gall formation in tomato plants.

Reference

1. Echeverrigaray, S., Zacaria, J., and Beltrão, R. Nematicidal activity of monoterpenoids against the root-knot nematode *Meloidogyne incognita*. *Phytopathology* **100**(2), 199-203 (2010).

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, 05/06/2021

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