

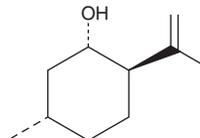
PRODUCT INFORMATION



(+)-Isopulegol

Item No. 23160

CAS Registry No.: 104870-56-6
Formal Name: 5S-methyl-2R-(1S-methylethenyl)-cyclohexanol
Synonym: d-Isopulegol
MF: C₁₀H₁₈O
FW: 154.3
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

(+)-Isopulegol is supplied as a liquid. A stock solution may be made by dissolving the (+)-isopulegol in water. The solubility of (+)-isopulegol in water is approximately 6 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(+)-Isopulegol is a terpenoid that has been found in *Cannabis*.¹

Reference

1. Fletcher, R.S. and McKay, J. Industrial hemp cannabis cultivars and seeds with stable cannabinoid profiles. *New West Genetics* **US2017/0339907A1** (2017).

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, 07/13/2018

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