

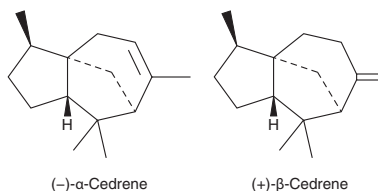
PRODUCT INFORMATION



Cedrene

Item No. 23180

CAS Registry No.: 11028-42-5
MF: $C_{15}H_{24}$
FW: 204.3
Purity: $\geq 95\%$ (mixture of isomers)
Supplied as: A neat oil
Storage: $4^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Cedrene is supplied as a neat oil. A stock solution may be made by dissolving the cedrene in the solvent of choice. Cedrene is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of cedrene in these solvents is approximately 20 mg/ml.

Cedrene is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cedrene should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Cedrene has a solubility of approximately 0.33 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cedrene is a sesquiterpene that has been found in a variety of plants, including *Cannabis*.¹

Reference

1. Elzinga, S., Fishedick, J., Podkolinski, R., *et al.* Cannabinoids and terpenes as chemotaxonomic markers in *Cannabis*. *Nat. Prod. Chem. Res.* **3**(4), 181 (2015).

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, 11/28/2022

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