

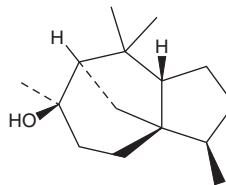
PRODUCT INFORMATION



(+)-Cedrol

Item No. 25769

CAS Registry No.: 77-53-2
Formal Name: (3R,3aS,6R,7R,8aS)-octahydro-3,6,8,8-tetramethyl-1H-3a,7-methanoazulen-6-ol
Synonym: α -Cedrol
MF: $C_{15}H_{26}O$
FW: 222.4
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

(+)-Cedrol is supplied as a solid. A stock solution may be made by dissolving the (+)-cedrol in the solvent of choice, which should be purged with an inert gas. (+)-Cedrol soluble in methanol (warmed) and is slightly soluble in chloroform.

Description

(+)-Cedrol is a sesquiterpene alcohol that has been used in *Cannabis* testing and has diverse biological activities.¹ It inhibits the growth of *L. sulphureus*, *G. trabeum*, *L. betulina*, and *T. versicolor* wood decay fungi when used at a concentration of 100 $\mu\text{g/ml}$.² (+)-Cedrol inhibits the cytochrome P450 (CYP) isoforms CYP2B6 and CYP3A4 ($K_s = 0.9$ and $3.4 \mu\text{M}$, respectively).³ *In vivo*, (+)-cedrol (200 mg/kg) prevents hair follicle dystrophy and reduces hair loss in a mouse model of alopecia induced by cyclophosphamide (Item No. 13849).⁴ (+)-Cedrol-loaded nanostructured lipid particles reduce peritoneal mast cell degranulation, a component of the type I anaphylactic response, induced by compound 48/80 (Item No. 22173) in mice.⁵

References

1. Fishedick, J.T. Identification of terpenoid chemotypes among high (-)-trans- Δ^9 -tetrahydrocannabinol-producing *Cannabis sativa* L. cultivars. *Cannabis Cannabinoid Res.* **2**(1), 34-47 (2017).
2. Cheng, S.S., Chung, M.J., Lin, C.Y., et al. Phytochemicals from *Cunninghamia konishii* Hayata act as antifungal agents. *J. Agric. Food Chem.* **60**(1), 124-128 (2012).
3. Jeong, H.U., Kwon, S.S., Kong, T.Y., et al. Inhibitory effects of cedrol, β -cedrene, and thujopsene on cytochrome P450 enzyme activities in human liver microsomes. *J. Toxicol. Environ. Health A.* **77**(22-24), 1522-1532 (2014).
4. Chen, S.S., Zhang, Y., Lu, Q.L., et al. Preventive effects of cedrol against alopecia in cyclophosphamide-treated mice. *Environ. Toxicol. Pharmacol.* **46**, 270-276 (2016).
5. Chakraborty, S., Kar, N., Kumari, L., et al. Inhibitory effect of a new orally active cedrol-loaded nanostructured lipid carrier on compound 48/80-induced mast cell degranulation and anaphylactic shock in mice. *Int. J. Nanomedicine* **12**, 4849-4868 (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, 09/29/2022

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