

PRODUCT INFORMATION



(-)-CP 47,497

Item No. 10917

CAS Registry No.: 114753-51-4
Formal Name: 2[(1S,3R)-3-hydroxycyclohexyl]-5-(2-methyloctan-2-yl)phenol

MF: C₂₁H₃₄O₂

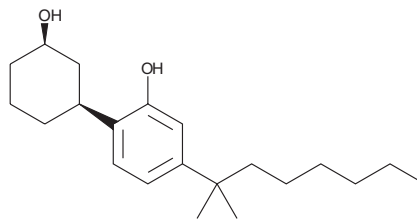
FW: 318.5

Purity: ≥98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Description

(-)-CP 47,497 (Item No. 10917) is an analytical reference standard that is structurally categorized as a cannabinoid (CB). It avidly binds the CB₁ receptor (K_i = 2.1 nM).¹ (-)-CP 47,497 is comparable or more potent than Δ⁹-THC (Item Nos. 12068 | ISO60157) in analgesic, motor depressant, anticonvulsant, and hypothermic effects in mice, rats, and dogs.^{2,3} (-)-CP 47,497 is regulated as a schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Melvin, L.S., Milne, G.M., Johnson, M.R., *et al.* Structure-activity relationships defining the ACD-tricyclic cannabinoids: Cannabinoid receptor binding and analgesic activity. *Drug Design Discov.* **13**, 155-166 (1995).
2. Melvin, L.S., Johnson, M.R., Harbert, C.A., *et al.* A cannabinoid derived prototypical analgesic. *J. Med. Chem.* **27**, 67-71 (1984).
3. Weissman, A., Milne, G.M., and Melvin, L.S., Jr. Cannabimimetic activity from CP-47,497, a derivative of 3-phenylcyclohexanol. *J. Pharmacol. Exp. Ther.* **223**(2), 516-523 (1982).

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, 03/06/2017

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