

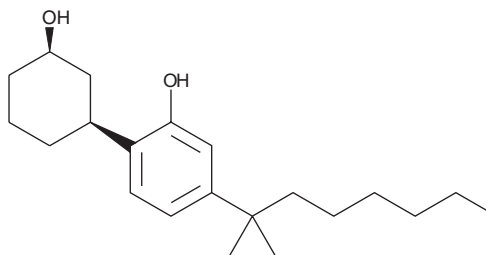
PRODUCT INFORMATION



(-)-CP 47,497 (exempt preparation)

Item No. 13218

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%
UV/Vis.: λ_{max}: 203, 218, 276 nm
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥3 years



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

Description

CP 47,497 is a bicyclic cannabinoid analog with potent analgesic activity.¹ It is comparable or more potent than Δ⁹-tetrahydrocannabinol in analgesic, motor depressant, anticonvulsant and hypothermic effects in mice, rats, and dogs.² The levorotatory enantiomer, (-)-CP 47,497 avidly binds the CB₁ receptor (K_i = 2.1 nM).³ (-)-CP 47,497 (exempt preparation) (Item No. 13218) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. Melvin, L.S., Johnson, M.R., Harbert, C.A., *et al.* A cannabinoid derived prototypical analgesic. *J. Med. Chem.* **27**(1), 67-71 (1984).
2. 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).
3. 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 Des. Discov.* **13**, 155-166 (1995).

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, 05/01/2026

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