

PRODUCT INFORMATION



(±)-CP 47,497

Item No. 10913

CAS Registry No.: 70434-82-1

Formal Name: *rel*-5-(1,1-dimethylheptyl)-2-[(1*R*,3*S*)-3-hydroxycyclohexyl]-phenol

Synonym: CP 47,947

MF: C₂₁H₃₄O₂

FW: 318.5

Purity: ≥95%

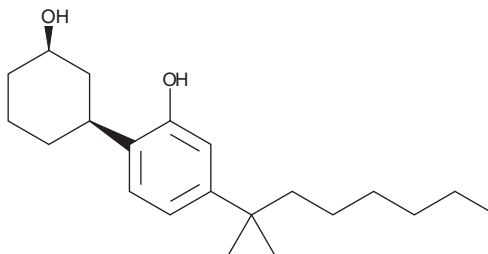
UV/Vis.: λ_{max}: 218, 276 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years

Packaged by: Cayman Chemical Company



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

Description

(±)-CP 47,497 is a monophenol cannabimimetic compound that binds the central cannabinoid (CB₁) receptor with a K_i value of 2.2 nM.^{1,2} It is equivalent in analgesic potency to D⁹-THC and exhibits other CB biological activities as well.

References

1. Huffman, J.W., Thompson, A.L.S., Wiley, J.L., *et al.* Synthesis and pharmacology of 1-deoxy analogs of CP-47,497 and CP-55,940. *Bioorg. Med. Chem.* **16**, 322-335 (2008).
2. Melvin, L.S., Milne, G.M., Johnson, M.R., *et al.* Structure-activity relationships for cannabinoid receptor-binding and analgesic activity: Studies of bicyclic cannabinoid analogs. *Mol. Pharmacol.* **44**, 1008-1015 (1993).

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, 07/13/2023

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