

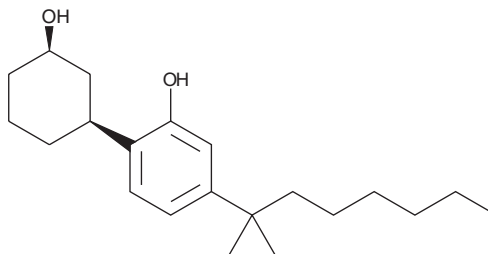
PRODUCT INFORMATION



(±)-CP 47,497 (exempt preparation)

Item No. 16851

CAS Registry No.: 70434-82-1
Formal Name: *rel*-2-[(1*S*,3*R*)-3-hydroxycyclohexyl]-5-(2-methyloctan-2-yl)-phenol
Synonym: CP 47,947
MF: C₂₁H₃₄O₂
FW: 318.5
Purity: ≥95%
UV/Vis.: λ_{max}: 218, 276 nm
Supplied as: A 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 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 Δ⁹-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**(1), 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**(5), 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, 01/15/2025

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