

# PRODUCT INFORMATION



## (±)-*epi* CP 47,497 (exempt preparation)

Item No. 13801

CAS Registry No.: 70434-83-2

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

Synonym: 3-*trans* CP 47,497

MF: C<sub>21</sub>H<sub>34</sub>O<sub>2</sub>

FW: 318.5

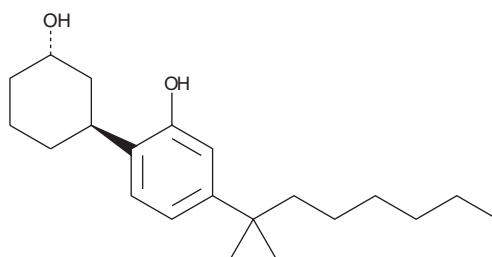
Purity: ≥98%

UV/Vis.: λ<sub>max</sub>: 219, 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

(±)-*epi* CP 47,497 (exempt preparation) (Item No. 13801) is an analytical reference standard that is structurally categorized as a cannabinoid (CB). It is an epimer of (±)-CP 47,497 (Item Nos. 10913 | 16851). (±)-CP 47,497 is a synthetic cannabinoid that avidly binds the CB<sub>1</sub> receptor (K<sub>i</sub> = 2.2 nM) and is equivalent in analgesic potency to Δ<sup>9</sup>-THC.<sup>1,2</sup> (±)-*epi* CP 47,497 is regulated as a Schedule 1 drug in the United States. (±)-*epi* CP 47,497 (exempt preparation) (Item No. 13801) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

### 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, 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