

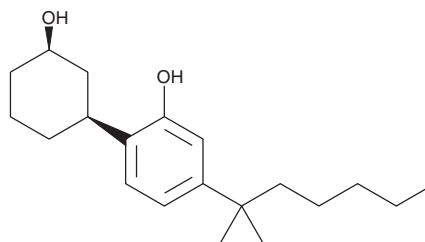
PRODUCT INFORMATION



CP 47,497-C6-homolog

Item No. 9001448

CAS Registry No.: 70435-06-2
Formal Name: cis-5-(1,1-dimethylhexyl)-2-(3-hydroxycyclohexyl)-phenol
MF: C₂₀H₃₂O₂
FW: 304.5
Purity: ≥95%
UV/Vis.: λ_{max}: 219, 276 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

CP 47,497-C6-homolog is a synthetic cannabinoid (CB) whose alkyl chain consists of six carbons. While *in vitro* data suggest that alkylaminoindoles with a side chain length of C4 to C6 possess equal or higher affinity to the CB central CB₁ and peripheral CB₂ receptors compared to Δ⁹-THC, the properties of this compound have not been evaluated.^{1,2} This product is intended for research and forensic applications.

References

1. Lindigkeit, R., Boehme, A., Eiserloh, I., *et al.* Spice: A never ending story? *Forensic Sci. Int.* **191(1-3)**, 58-63 (2009).
2. 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).

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, 07/11/2023

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