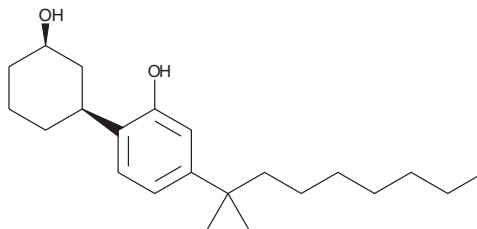


PRODUCT INFORMATION



(±)-CP 47,497-C8-homolog Item No. 10908

CAS Registry No.: 70434-92-3
Formal Name: *rel*-5-(1,1-dimethyloctyl)-2-[(1*R*,3*S*)-3-hydroxycyclohexyl]-phenol
Synonyms: Cannabicyclohexanol, CAY10596
MF: C₂₂H₃₆O₂
FW: 332.5
Purity: ≥98%
UV/Vis.: λ_{max}: 277 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Description

(±)-CP 47,497-C8-homolog (Item No. 10908) is an analytical reference standard categorized as a synthetic cannabinoid.^{1,2} CP 47,497-C8-homolog has been found in Spice herbal blends.³ (±)-CP 47,497-C8-homolog is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. 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).
2. Compton, D.R., Johnson, M.R., Melvin, L.S., *et al.* Pharmacological profile of a series of bicyclic cannabinoid analogs: Classification as cannabimimetic agents. *J. Pharmacol. Exp. Ther.* **260**(1), 201-209 (1992).
3. Gwak, S., Arroyo-Mora, L.E., and Almirall, J.R. Qualitative analysis of seized synthetic cannabinoids and synthetic cathinones by gas chromatography triple quadrupole tandem mass spectrometry. *Drug Test Anal.* **7**(2), 131-130 (2015).

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, 01/08/2018

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