

PRODUCT INFORMATION



(±)-9β-hydroxy Hexahydrocannabinol

Item No. 35266

Formal Name: (±)-rel-(6aR,9R,10aR)-6,6,9-trimethyl-3-pentyl-6a,7,8,9,10,10a-hexahydro-6H-benzo[c]chromene-1,9-diol

Synonym: (±)-9β-hydroxy HHC

MF: C₂₁H₃₂O₃

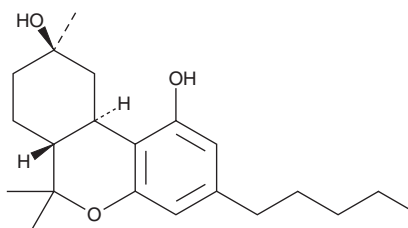
FW: 332.5

Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



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

Description

(±)-9β-hydroxy Hexahydrocannabinol (Item No. 35266) is an analytical reference standard that is structurally similar to known phytocannabinoids.¹ (±)-9β-hydroxy Hexahydrocannabinol is a metabolite of 9-hexahydrocannabinol.² It has been found as an impurity in products marketed as containing Δ⁸-THC.³ This product is intended for research and forensic applications.

Reference

1. Nalli, Y., Jan, S., Lauro, G., *et al.* Isolation, synthesis and structure determination of cannabidiol derivatives and their cytotoxic activities. *Nat. Prod. Res.* **35**(3), 471-480 (2021).
2. Kobidze, G., Sprega, G., Montanari, E., *et al.* The first LC-MS/MS stereoselective bioanalytical methods to quantitatively detect 9R- and 9S-hexahydrocannabinols and their metabolites in human blood, oral fluid and urine. *J. Pharm. Biomed. Anal.* **240**, 115918 (2024).
3. Radwan, M.M., Wanas, A.S., Gul, W., *et al.* Isolation and characterization of impurities in commercially marketed Δ⁸-THC products. *J. Nat. Prod.* **86**(4), 822-829 (2023).

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, 03/15/2024

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