

PRODUCT INFORMATION



(±)-9α-hydroxy Hexahydrocannabinol

Item No. 36129

CAS Registry No.: 16964-49-1

Formal Name: (6aR,9S,10aR)-*rel*-6a,7,8,9,10,10a-hexahydro-6,6,9-trimethyl-3-pentyl-6H-dibenzo[b,d]pyran-1,9-diol

Synonyms: (±)-9α-hydroxy HHC, (±)-9α-OH-HHC

MF: C₂₁H₃₂O₃

FW: 332.5

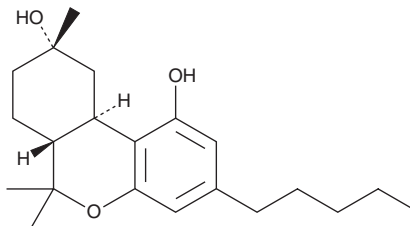
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥5 years

Item Origin: Synthetic



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

Description

(±)-9α-hydroxy Hexahydrocannabinol (Item No. 36129) is an analytical reference standard that is structurally similar to known phytocannabinoids.¹ (±)-9α-hydroxy Hexahydrocannabinol is a metabolite of 9-hexahydrocannabinol.² It has been found in high potency strains of *C. sativa*.¹ (±)-9α-hydroxy Hexahydrocannabinol has also been found as an impurity in products marketed as containing Δ⁸-THC.³ (±)-9α-hydroxy Hexahydrocannabinol is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Ahmed, S.A., Ross, S.A., Slade, D., *et al.* Minor oxygenated cannabinoids from high potency *Cannabis sativa* L. *Phytochemistry* **117**, 194-199 (2015).
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**, 15918 (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, 06/25/2025

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