

PRODUCT INFORMATION



11-hydroxy-9(R)-Hexahydrocannabinol

Item No. 36353

CAS Registry No.: 64663-39-4

Formal Name: 6aR,7,8,9R,10,10aR-hexahydro-1-hydroxy-6,6-dimethyl-3-pentyl-6H-dibenzo[b,d]pyran-9-methanol

Synonyms: 11-hydroxy-9 β -Hexahydrocannabinol, 11-hydroxy-9 β -HHC, 11-hydroxy-9(R)-HHC

MF: C₂₁H₃₂O₃

FW: 332.5

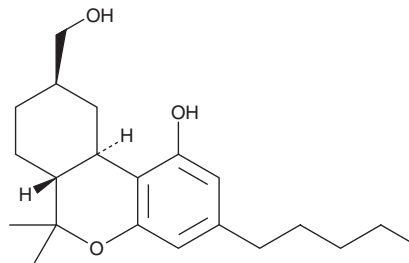
Purity: $\geq 95\%$

Supplied as: A neat solid

Storage: -20°C

Stability: ≥ 5 years

Item Origin: Plant/Semi-synthetic



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

Description

11-hydroxy-9(R)-Hexahydrocannabinol (Item No. 36353) is an analytical reference standard categorized as a phytocannabinoid metabolite.¹⁻⁴ 11-hydroxy-9(R)-Hexahydrocannabinol is an active metabolite of exo-THC (Item No. 36898), 9-hexahydrocannabinol, and Δ^9 -THC (Item Nos. ISO60157 | 9003771 | 12068).¹⁻³ It induces generalization to Δ^9 -THC in drug discrimination tests in rats and pigeons.⁴ 11-hydroxy-9(R)-Hexahydrocannabinol is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Harvey, D.J., Gill, E.E., Slater, M., *et al.* Identification of the *in vivo* liver metabolites of (-)- Δ^7 -tetrahydrocannabinol produced by the mouse. *Drug Metab. Dispos.* **8**(6), 439-445 (1980).
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. Jørgensen, C.F., Rasmussen, B.S., Linnet, K., *et al.* Evidence of 11-hydroxy-hexahydrocannabinol and 11-nor-9-carboxy-hexahydrocannabinol as Novel Human Metabolites of Δ^9 -tetrahydrocannabinol. *Metabolites* **13**(12), 1169 (2023).
4. Järbe, T.U., Hiltunen, A.J., Lander, N., *et al.* Cannabimimetic activity (Δ^1 -THC cue) of cannabidiol monomethyl ether and two stereoisomeric hexahydrocannabinols in rats and pigeons. *Pharmacol. Biochem. Behav.* **25**(2), 393-399 (1986).

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