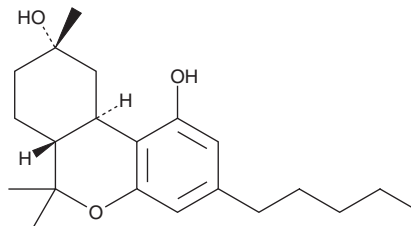


PRODUCT INFORMATION

(±)-9α-hydroxy Hexahydrocannabinol (exempt preparation)

Item No. 44126

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
Synonym: (±)-9α-hydroxy HHC
MF: C₂₁H₃₂O₃
FW: 332.5
Purity: ≥98%
Supplied as: A 1 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥3 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. 44126) 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.¹ (±)-9α-hydroxy Hexahydrocannabinol is regulated as a Schedule I compound in the United States. (±)-9α-hydroxy Hexahydrocannabinol (Item No. 44126) is provided as a DEA exempt preparation. 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. 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, 06/27/2025

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