

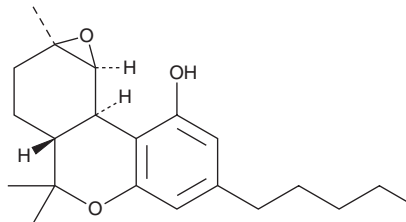
PRODUCT INFORMATION



9 β ,10 β -epoxy Hexahydrocannabinol

Item No. 42023

CAS Registry No.: 1781270-07-2
Formal Name: 1aR,2,3aR,4,9bR,9cS-hexahydro-1a,4,4-trimethyl-7-pentyl-3H-oxireno[3,4]benzo[1,2-c][1]benzopyran-9-ol
Synonyms: 9 β ,10 β -EHHHC, 9 β ,10 β -epoxy HHC
MF: C₂₁H₃₀O₃
FW: 330.5
Purity: $\geq$ 95%
Supplied as: A 1 mg/ml solution in acetonitrile
Storage: -20°C
Stability: $\geq$ 2 years



Special Conditions: Protect from light, moisture, and oxygen

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

Description

9 β ,10 β -epoxy Hexahydrocannabinol (Item No. 42023) is an analytical reference standard categorized as a phytocannabinoid.¹ 9 β ,10 β -epoxy Hexahydrocannabinol induces catalepsy, hypothermia, hypolocomotion, and antinociception in the mouse tetrad assay for cannabimimetic activity. It is a potential metabolite of Δ^9 -THC (Item Nos. ISO60157 | 42138 | 12068) based on the published metabolism of Δ^8 -THC (Item Nos. ISO60158 | 41193 | 14042).² At the time 9 β ,10 β -epoxy hexahydrocannabinol (Item No. 42023) was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. 9 β ,10 β -epoxy Hexahydrocannabinol is regulated as a Schedule I product. This product is intended for research and forensic applications.

References

1. Radwan, M.M., ElSohly, M.A., El-Alfry, A.T., *et al.* Isolation and pharmacological evaluation of minor cannabinoids from high-potency Cannabis sativa. *J. Nat. Prod.* **78**(6), 1271-1276 (2015).
2. Narimatsu, S., Watanabe, K., Shimonishi, T., *et al.* Metabolic disposition of 8 α , 9 α - and 8 β , 9 β -epoxyhexahydrocannabinols in the mouse. *J. Pharmacobiodyn.* **7**(9), 671-676 (1984).

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, 11/24/2025

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