

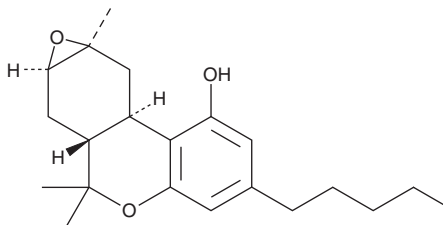
PRODUCT INFORMATION



8β,9β-epoxy Hexahydrocannabinol (exempt preparation)

Item No. 44608

CAS Registry No.: 64783-83-1
Formal Name: 6aR,7,7aR,8aS,9,9aR-hexahydro-6,6,8a-trimethyl-3-pentyl-6H-oxireno[4,5]benzo[1,2-c][1]benzo-1-ol
Synonyms: 8β,9β-EHHC, 8β,9β-epoxy HHC
MF: C₂₁H₃₀O₃
FW: 330.5
Purity: ≥95%
Supplied as: A 1 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥3 years



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

Description

8β,9β-epoxy Hexahydrocannabinol (Item No. 44608) is an analytical reference standard categorized as a phytocannabinoid metabolite.¹ 8β,9β-epoxy Hexahydrocannabinol is an active metabolite of Δ⁸-THC (Item Nos. ISO60158 | 41193 | 14042). It prolongs pentobarbital-induced sleeping time in mice. 8β,9β-epoxy Hexahydrocannabinol is regulated as a Schedule I compound in the United States. 8β,9β-epoxy Hexahydrocannabinol (exempt preparation) (Item No. 44608) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

Reference

1. Narimatsu, S., Watanabe, K., Shimonishi, T., *et al.* Metabolic disposition of 8α, 9α- and 8β, 9β-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, 09/19/2025

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