

PRODUCT INFORMATION

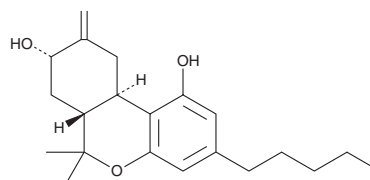


8 α -hydroxy-exo-THC (exempt preparation)

Item No. 39795

Formal Name: 6,6-dimethyl-9-methylene-3-pentyl-6aR,7,8S,9,10,10aR-hexahydro-6H-benzo[c]chromene-1,8-diol

Synonyms: 6 α -hydroxy- $\Delta^{1(7)}$ -Tetrahydrocannabinol, 6(S)-hydroxy- $\Delta^{1(7)}$ -Tetrahydrocannabinol, 6 α -hydroxy- $\Delta^{1(7)}$ -THC, 6(S)-hydroxy- $\Delta^{1(7)}$ -THC, 8 α -hydroxy- $\Delta^{9(11)}$ -Tetrahydrocannabinol, 8(S)-hydroxy- $\Delta^{9(11)}$ -Tetrahydrocannabinol, 8 α -hydroxy- $\Delta^{9(11)}$ -THC, 8(S)-hydroxy- $\Delta^{9(11)}$ -THC, 8 α -hydroxy-exo-Tetrahydrocannabinol, 8(S)-hydroxy-exo-Tetrahydrocannabinol, 8 α -hydroxy-exo-THC, 8(S)-hydroxy-exo-THC



MF: C₂₁H₃₀O₃
FW: 330.5
Purity: ≥98%
Supplied as: A 100 µg/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

8 α -hydroxy-exo-THC (Item No. 39795) (exempt preparation) is an analytical reference standard that is structurally similar to known phytocannabinoids. 8 α -hydroxy-exo-THC is a metabolite of exo-THC (Item No. 36898).¹ 8 α -hydroxy-exo-THC is regulated as a Schedule I compound in the United States. 8 α -hydroxy-exo-THC (exempt preparation) (Item No. 39795) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

Reference

1. Harvey, D.J. and Brown, N.K. Comparative in vitro metabolism of the cannabinoids. *Pharmacol. Biochem. Behav.* **40**(3), 533-540 (1991).

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, 01/15/2025

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