

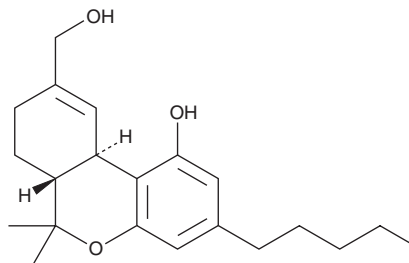
PRODUCT INFORMATION



(-)-11-hydroxy- Δ^9 -THC

Item No. 14239

CAS Registry No.: 36557-05-8
Formal Name: 6aR,7,8,10aR-tetrahydro-1-hydroxy-6,6-dimethyl-3-pentyl-6H-dibenzo[b,d]pyran-9-methanol
Synonyms: (-)-7-hydroxy- Δ^1 -Tetrahydrocannabinol, (-)-11-hydroxy- Δ^9 -Tetrahydrocannabinol, (-)-7-hydroxy- Δ^1 -THC
MF: C₂₁H₃₀O₃
FW: 330.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Description

(-)-11-hydroxy- Δ^9 -THC (Item No. 14239) is an analytical reference standard categorized as a phytocannabinoid metabolite.¹⁻⁴ (-)-11-hydroxy- Δ^9 -THC is a metabolite of Δ^9 -THC (Item Nos. ISO60157 | 9003771 | 12068). (-)-11-hydroxy- Δ^9 -THC is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Martin, B.R., Balster, R.L., Razdan, R.K., *et al.* Behavioral comparisons of the stereoisomers of tetrahydrocannabinols. *Life Sci.* **29(6)**, 565-574 (1981).
2. Grafinger, K.E. and Weinmann, W. Determination of the cross-reactivity of the biological metabolite (-)-*trans*- Δ^9 -tetrahydrocannabinol-carboxylic acid-glucuronide (THC-COOH-Gluc) for cannabinoid immunoassays. *J. Anal. Toxicol.* **45(3)**, 291-296 (2021).
3. Bland, T.M., Haining, R.L., Tracy, T.S., *et al.* CYP2C-catalyzed delta(9)-tetrahydrocannabinol metabolism: Kinetics, pharmacogenetics and interaction with phenytoin. *Biochem. Pharmacol.* **70(7)**, 1096-1103 (2005).
4. Coles, R., Clements, T.T., Nelson, G.J., *et al.* Simultaneous analysis of the Δ^9 -THC metabolites 11-nor-9-carboxy- Δ^9 -THC and 11-hydroxy- Δ^9 -THC in meconium by GC-MS. *J. Anal. Toxicol.* **29(6)**, 522-527 (2016).

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, 04/24/2024

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