

PRODUCT INFORMATION



(-)-11-hydroxy- Δ^8 -Tetrahydrocannabivarin

Item No. 36630

CAS Registry No.: 1421594-84-4
Formal Name: 6aR,7,10,10aR-tetrahydro-1-hydroxy-6,6-dimethyl-3-propyl-6H-dibenzo[b,d]pyran-9-methanol

Synonyms: (-)-11-hydroxy- Δ^8 -THCV,
(-)-11-hydroxy- Δ^8 -THV

MF: C₁₉H₂₆O₃

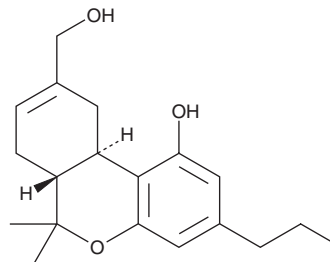
FW: 302.4

Purity: \geq 98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: \geq 4 years



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

Description

(-)-11-hydroxy- Δ^8 -Tetrahydrocannabivarin (Item No. 36630) is an analytical reference standard categorized as a synthetic cannabinoid.¹ (-)-11-hydroxy- Δ^8 -Tetrahydrocannabivarin is a metabolite of Δ^8 -tetrahydrocannabivarin (Item Nos. 27322 | 38795). (-)-11-hydroxy- Δ^8 -Tetrahydrocannabivarin is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

Reference

1. Bátkai, S., Mukhopadhyay, P., Horváth, B., *et al.* Δ^8 -Tetrahydrocannabivarin prevents hepatic ischaemia/reperfusion injury by decreasing oxidative stress and inflammatory responses through cannabinoid CB₂ receptors. *Br. J. Pharmacol.* **165**(8), 2450-2461 (2012).

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, 07/19/2023

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