

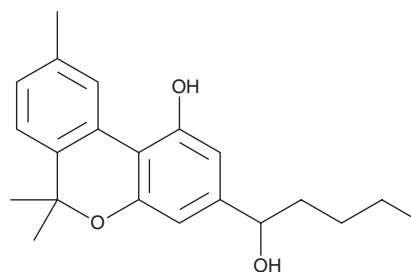
PRODUCT INFORMATION



(±)-1'-hydroxy Cannabinol

Item No. 44316

CAS Registry No.: 64663-47-4
Formal Name: α-butyl-1-hydroxy-6,6,9-trimethyl-6H-dibenzo[b,d]pyran-3-methanol
Synonyms: CBN-1'-hydroxy, 1'-hydroxy CBN
MF: C₂₁H₂₆O₃
FW: 326.4
Purity: ≥98%
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

(±)-1'-hydroxy Cannabinol (Item No. 44316) is an analytical reference standard categorized as a phytocannabinoid metabolite.¹ (±)-1'-hydroxy Cannabinol is a minor metabolite of cannabinol (Item Nos. ISO60183 | 42139 | 25495). This product is intended for research and forensic applications.

Reference

1. Kundu, D., Franchini, L., Hasdemir, H.S., *et al.* Distinct interactions of cannabinol and its cytochrome P450-generated metabolites with receptors and sensory neurons. *J. Med. Chem.* **68(13)**, 13935-13953 (2025).

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, 09/03/2025

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