

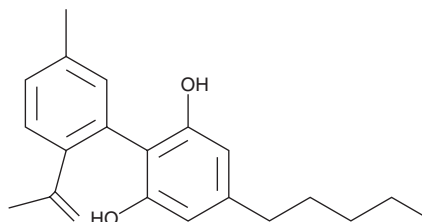
PRODUCT INFORMATION



Cannabinodiol

Item No. 9002479

CAS Registry No.: 39624-81-2
Formal Name: 5'-methyl-2'-(1-methylethenyl)-4-pentyl-[1,1'-biphenyl]-2,6-diol
Synonym: CBND
MF: C₂₁H₂₆O₂
FW: 310.4
Purity: ≥98%
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Description

Cannabinodiol (Item No. 9002479) is an analytical reference standard that is categorized as a phytocannabinoid. It is an aromatic compound that has been isolated from Lebanese hashish.¹ The physiological and toxicological properties of this compound are not known. This product is intended for research and forensic applications.

Reference

1. Lousberg, R.J.J., Bercht, C.A.L., van Ooyen, R., *et al.* Cannabinodiol: Conclusive identification and synthesis of a new cannabinoid from *Cannabis sativa*. *Phytochem.* **16(5)**, 595-597 (1977).

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, 07/25/2023

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