

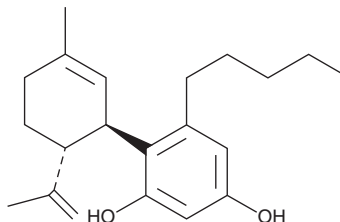
PRODUCT INFORMATION



Abnormal Cannabidiol

Item No. 10004259

CAS Registry No.: 22972-55-0
Formal Name: 4-[(1R,6R)-3-methyl-6-(1-methylethenyl)-2-cyclohexen-1-yl]-5-pentyl-1,3-benzenediol
Synonyms: Abn-CBD, *ortho*-Cannabidiol, *o*-CBD
MF: C₂₁H₃₀O₂
FW: 314.5
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥4 years



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

Description

Abnormal cannabidiol (Item No. 10004259) is an analytical reference standard that is structurally similar to known phytocannabinoids. It is a bioproduct in the synthesis of cannabidiol.¹ Abnormal cannabidiol is not found in Cannabis and has been used as an internal standard for the detection of cannabinoids in hemp extracts.² This product is intended for research and forensic applications.

References

1. Chiurchiù, E., Sampaolesi, S., Allegrini, P., *et al.* A novel and practical continuous flow chemical synthesis of cannabidiol (CBD) and its CBDV and CBDB analogues. *Eur. J. Org. Chem.* **8**, 1286-1289 (2021).
2. Song, L., Carlson, S., Valenzuela, G., *et al.* Development of a validated method for rapid quantification of up to sixteen cannabinoids using ultra-high-performance liquid chromatography diode-array detector with optional electrospray ionization time-of-flight mass spectrometry detection. *J. Chromatogr. A* **1670**, 462953 (2022).

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, 07/07/2023

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