

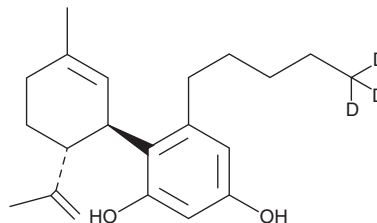
PRODUCT INFORMATION



Abnormal Cannabidiol-d₃

Item No. 10009523

CAS Registry No.: 2649096-15-9
Formal Name: 4-[(1R,6R)-3-methyl-6-(1-methylethenyl)-2-cyclohexen-1-yl]-5-pentyl-5,5,5-d₃-1,3-benzenediol
Synonyms: Abn-CBD-d₃, o-CBD, *ortho*-Cannabidiol
MF: C₂₁H₂₇D₃O₂
FW: 317.5
Chemical Purity: ≥98% (Abnormal cannabidiol)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A 1 mg/ml 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-d₃ (Item No. 10009523) is intended for use as an internal standard for the quantification of abnormal cannabidiol (Item No. 10004259) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Abnormal cannabidiol 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, 10/27/2025

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