

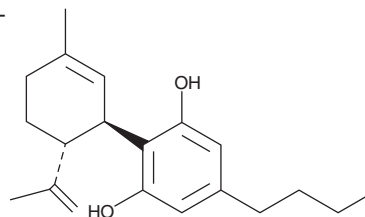
PRODUCT INFORMATION



CBDB

Item No. 29031

CAS Registry No.: 60113-11-3
Formal Name: (1*R-trans*)-5-butyl-2-[3-methyl-6-(1-methylethenyl)-2-cyclohexen-1-yl]-1,3-benzenediol
Synonyms: Cannabidibutol, Cannabidiol-C4, CBD-C4
MF: C₂₀H₂₈O₂
FW: 300.4
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

CBDB (Item No. 29031) is an analytical reference standard categorized as a phytocannabinoid.¹ CBDB has been found in strains of *Cannabis* and hemp.^{1,2} CBDB is considered an impurity in commercial extractions of cannabidiol (CBD) from hemp.² This product is intended for research and forensic application

References

1. Linciano, P., Citti, C., Luongo, L., *et al.* Isolation of a high-affinity cannabinoid for the human CB₁ receptor from a medicinal *Cannabis sativa* variety: Δ⁹-Tetrahydrocannabutol, the butyl homologue of Δ⁹-tetrahydrocannabinol. *J. Nat. Prod.* **83**(1), 88-98 (2019).
2. Citti, C., Linciano, P., Forni, F., *et al.* Analysis of impurities of cannabidiol from hemp. Isolation, characterization and synthesis of *cannabidibutol*, the novel cannabidiol butyl analog. *J. Pharm. Biomed. Anal.* **175**, 112752 (2019).

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, 01/14/2026

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