

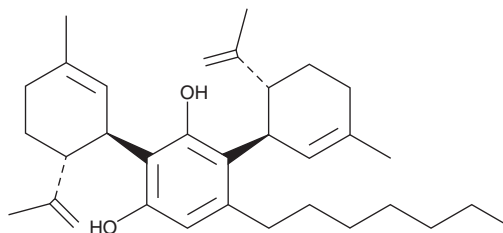
PRODUCT INFORMATION



di-coupled CBDP

Item No. 41323

CAS Registry No.: 1260669-31-5
Formal Name: 5-heptyl-2,4-bis[(1R)-3-methyl-6R-(1-methylethenyl)-2-cyclohexen-1-yl]-1,3-benzenediol
Synonyms: bis-menthadiene Cannabidiphorol, di-coupled Cannabidiphorol, di-coupled CBD-C7, di-coupled CBD-heptyl, bis-CBDP isomer, dialkylated CBDP
MF: C₃₃H₄₈O₂
FW: 476.7
Purity: ≥95%
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

di-coupled CBDP (Item No. 41323) is an analytical reference standard that is structurally similar to known phytocannabinoids. di-coupled CBDP is a dialkylated form of CBDP (Item Nos. 33611 | 30169) and a byproduct formed in the synthesis of CBDP.¹ This product is intended for research and forensic applications.

Reference

1. Golliher, A.E., Tenorio, A.J., Dimauro, N.O., *et al.* Using (+)-carvone to access novel derivatives of (+)- *ent*-cannabidiol: The first asymmetric syntheses of (+)- *ent*-CBDP and (+)- *ent*-CBDV. *Tetrahedron Lett.* **67**, 152891 (2021).

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, 11/21/2025

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