

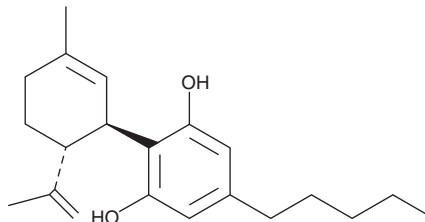
Product Information



Cannabidiol (DEA Schedule I Regulated Compound)

Catalog No. 90080 • Lot No. XXXXXX • Size: XX mg

NDC: 65326-002-XX
CAS Registry No.: 13956-29-1
Formal Name: 2-[1R-3-methyl-6R-(1-methylethenyl)-2-cyclohexen-1-yl]-5-pentyl-1,3-benzenediol
MF: C₂₁H₃₀O₂
FW: 314.5
Purity: ≥99%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid
Packaged by: Cayman Chemical Company



Laboratory Procedures

Cannabidiol is one of the major phenolic components of hemp. In contrast to the minor component Δ^9 -THC, it is not psychoactive. For long term storage, we suggest that cannabidiol be stored as supplied at -20°C. It will be stable for at least one year.

Cannabidiol is supplied as a crystalline solid. A stock solution may be made by dissolving the cannabidiol in an organic solvent purged with an inert gas. Cannabidiol is soluble in organic solvents such as ethanol (36 mg/ml), DMSO (60 mg/ml), and dimethyl formamide (50 mg/ml). It will be stable for at least six months in these solvents if stored at -20°C.

Cannabidiol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cannabidiol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Cannabidiol has a solubility of 250 µg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. Store aqueous solutions of cannabidiol on ice and use within 12 hours of preparation. Although the aqueous solutions of cannabidiol may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

Antioxidant activity has been reported to be a general property of the phenolic components of marijuana. Unlike Δ^9 -THC, cannabidiol can be administered at relatively high doses without undesired toxic or psychological effects.¹ Cannabidiol at a concentration of 10 µM was neuroprotective against both excitatory neurotransmitter (glutamate) and oxidant (hydroperoxide) induced neurotoxicity.²

References

1. Consroe, P., Laguna, J., Allender, J., *et al.* Controlled clinical trial of cannabidiol in Huntington's disease. *Pharmacol. Biochem. Behav.* **40**, 701-708 (1991).
2. Hampson, A.J., Grimaldi, M., Axelrod, J., *et al.* Cannabidiol and (-) Δ^9 -tetrahydrocannabinol are neuroprotective antioxidants. *Proc. Nat. Acad. Sci USA* **95**, 8268-8273 (1998).

Related Products

2-Arachidonoyl Glycerol - Cat. No. 62160 • Arachidonoyl Ethanolamide - Cat. No. 90050 • HU-210 - Cat. No. 90082 • CB₁ Receptor Polyclonal Antibody - Cat. No. 101500 • CB₂ Receptor Polyclonal Antibody - Cat. No. 101550

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 03/10/2009