

PRODUCT INFORMATION



Cannabigerolic Acid-d₉

Item No. 44742

Formal Name: 3-[(2E)-3,7-dimethyl-2,6-octadien-1-yl]-2,4-dihydroxy-6-(pentyl-2,2,3,3,4,4,5,5,5-d₉)benzoic acid

Synonym: CBGA-d₉

MF: C₂₂H₂₃D₉O₄

FW: 369.5

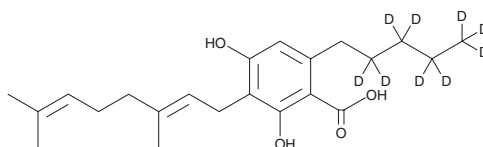
Chemical Purity: ≥98% (cannabigerolic acid)

Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀

Supplied as: A 1 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

Cannabigerolic acid-d₉ (Item No. 44742) is an analytical reference standard intended for use as an internal standard for the quantification of cannabigerolic acid (Item Nos. 20019 | 42523 | 9001572) 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).

Cannabigerolic acid is categorized as a phytocannabinoid.¹ It is a precursor in the biosynthesis of tetrahydrocannabinolic acid, cannabidiolic acid (Item Nos. 18090 | 41378 | 36871), and cannabichromenic acid (Item Nos. 37558 | 43097 | 36998). This product is intended for research and forensic applications.

Reference

1. Morimoto, S., Taura, F., and Shoyama, Y. Biosynthesis of cannabinoids in *Cannabis sativa* L. *Curr. Top. Phytochem.* 2, 103-113 (1999).

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, 06/18/2026

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