

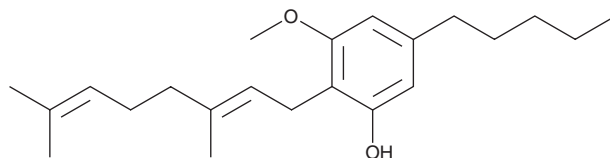
PRODUCT INFORMATION



CBGM

Item No. 9003893

CAS Registry No.: 29106-17-0
Formal Name: 2-[(2E)-3,7-dimethyl-2,6-octadien-1-yl]-3-methoxy-5-pentyl-phenol
Synonyms: O-methyl Cannabigerol, Cannabigerol monomethyl ether
MF: C₂₂H₃₄O₂
FW: 330.5
Purity: ≥98%
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Description

CBGM (Item No. 9003893) is an analytical reference standard categorized as a phytocannabinoid.¹ This product is intended for research and forensic applications.

Reference

1. Linciano, P., Citti, C., Russo, F., *et al.* Identification of a new cannabidiol n-hexyl homolog in a medicinal cannabis variety with an antinociceptive activity in mice: Cannabidihexol. *Sci. Rep.* **10(1)**, 22019 (2020).

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, 01/13/2026

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