

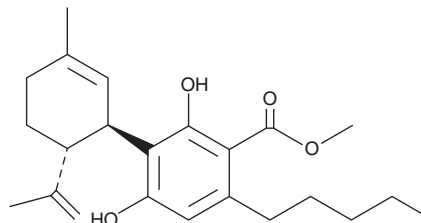
PRODUCT INFORMATION



Cannabidiolic Acid methyl ester

Item No. 28347

CAS Registry No.: 55658-71-4
Formal Name: 2,4-dihydroxy-3-[(1R,6R)-3-methyl-6-(1-methylethenyl)-2-cyclohexen-1-yl]-6-pentyl-benzoic acid, methyl ester
Synonyms: CBDA methyl ester, HU-580
MF: C₂₃H₃₂O₄
FW: 372.5
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 268 nm
Supplied as: A 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

Cannabidiolic acid methyl ester (Item No. 28347) is an analytical reference standard that is the methyl ester form of cannabidiolic acid (Item No. 18090). Cannabidiolic acid methyl ester has analgesic activity in rats.¹ This product is intended for research and forensic applications.

Reference

1. Zhu, Y.F., Linher-Melville, K., Niazmand, M.J., *et al.* An evaluation of the anti-hyperalgesic effects of cannabidiolic acid-methyl ester in a preclinical model of peripheral neuropathic pain. *Br. J. Pharmacol.* **177(12)**, 2712-2725 (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, 07/10/2023

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