

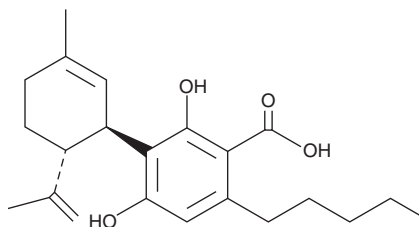
# PRODUCT INFORMATION



## Cannabidiolic Acid

Item No. 36871

**CAS Registry No.:** 1244-58-2  
**Formal Name:** 2,4-dihydroxy-3-[(1R,6R)-3-methyl-6-(1-methylethenyl)-2-cyclohexen-1-yl]-6-pentyl-benzoic acid  
**Synonyms:** CBDA  
**MF:** C<sub>22</sub>H<sub>30</sub>O<sub>4</sub>  
**FW:** 358.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 224, 270 nm  
**Supplied as:** A solid  
**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 (Item No. 36871) is an analytical reference standard categorized as a phytocannabinoid.<sup>1,2</sup> Cannabidiolic acid has anxiolytic-like and antinociceptive activities in mice.<sup>2,3</sup> This product is intended for research and forensic applications.

### References

1. Taura, F., Morimoto, S., and Shoyama, Y. Purification and characterization of cannabidiolic-acid synthase from *Cannabis sativa* L.. Biochemical analysis of a novel enzyme that catalyzes the oxidocyclization of cannabigerolic acid to cannabidiolic acid. *J. Biol. Chem.* **271(29)**, 17411-17416 (1996).
2. ZagZoog, A., Mohamed, K.A., Kim, H.J.J., et al. In vitro and in vivo pharmacological activity of minor cannabinoids isolated from *Cannabis sativa*. *Sci. Rep.* **10(1)**, 20405 (2020).
3. Vigli, D., Cosentino, L., Pellas, M., et al. Chronic treatment with cannabidiolic acid (CBDA) reduces thermal pain sensitivity in male mice and rescues the hyperalgesia in a mouse model of Rett syndrome. *Neuroscience* **453**, 113-123 (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, 01/08/2024

#### 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