

# PRODUCT INFORMATION



## $\Delta^9$ -THCB Butanoate

Item No. 43582

**Formal Name:** 3-butyl-6,6,9-trimethyl-6aR,7,8,10aR-tetrahydro-6H-benzo[c]chromen-1-yl butyrate

**Synonyms:**  $\Delta^9$ -Tetrahydrocannabinol-C4 Butanoate,  $\Delta^9$ -Tetrahydrocannabitol Butanoate,  $\Delta^9$ -THCB-O-Butanoate,  $\Delta^9$ -THCB Butyrate,  $\Delta^9$ -THC-C4 Butanoate

**MF:** C<sub>24</sub>H<sub>34</sub>O<sub>3</sub>

**FW:** 370.5

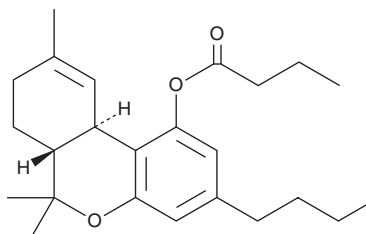
**Purity:**  $\geq 95\%$

**Supplied as:** A 1 mg/ml solution in 99:1 acetonitrile:DIPEA with 0.05% ascorbic acid

**Storage:** -80°C

**Stability:**  $\geq 2$  years

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



### Description

$\Delta^9$ -THCB butanoate (Item No. 43582) is an analytical reference standard that is structurally similar to known phytocannabinoids.  $\Delta^9$ -THCB butanoate has been found in electronic cigarette (e-cigarette) cartridge oil.<sup>1,2</sup>  $\Delta^9$ -THCB butanoate is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

### References

1. Tanaka, R., Ito, M., and Kikura-Hanajiri, R. Identification of acetyl, propionyl, and butanoyl derivatives of THC and its analogs, and their synthetic by-products in oils and a herbal product. *Drug Test. Anal.* **17(9)**, 1795-1802 (2025).
2. Tanaka, R., Kawamura, M., Ito, M., *et al.* Identification of novel THC analogs: Chloro derivatives of hexahydrocannabinol and tetrahydrocannabitol butanoate in an oil product obtained in Japan. *Drug Test Anal.* (2025).

#### 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, 12/15/2025

#### 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