

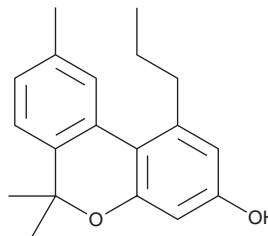
PRODUCT INFORMATION



Abnormal Cannabivarin

Item No. 36919

Formal Name: 6,6,9-trimethyl-1-propyl-6H-benzo[c]chromen-3-ol
Synonyms: Abn-CBV, Abnormal Cannabivarol
MF: C₁₉H₂₂O₂
FW: 282.4
Purity: ≥98%
UV/Vis.: λ_{max}: 280 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

Abnormal cannabivarin (Item No. 36919) is an analytical reference standard that is structurally similar to known phytocannabinoids. Abnormal cannabivarin is a potential regioisomer produced in the synthesis of cannabivarin (Item No. 21664) based on the published synthesis of Δ^8 -THC (Item No. 14042).¹ This product is intended for research and forensic applications.

Reference

1. Razdan, R.K., Dalzell, H.C., and Handrick, G.R. Hashish. A simple one-step synthesis of (-)- Δ^1 -tetrahydrocannabinol (THC) from p-mentha-2,8-dien-1-ol and olivetol. *J. Am. Chem. Soc.* **96(18)**, 5860-5865 (1974).

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, 07/19/2023

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