

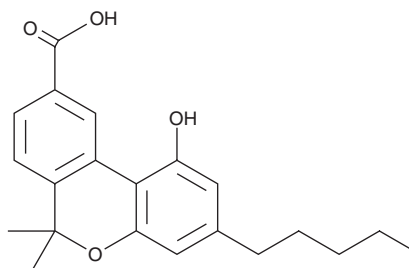
PRODUCT INFORMATION



11-nor-9-carboxy Cannabinol

Item No. 34592

CAS Registry No.: 53989-32-5
Formal Name: 1-hydroxy-6,6-dimethyl-3-pentyl-6H-dibenzo[b,d]pyran-9-carboxylic acid
Synonym: Cannabinol-11-oic Acid
MF: C₂₁H₂₄O₄
FW: 340.4
Purity: ≥95%
UV/Vis.: λ_{max}: 230, 287 nm
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥2 years



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

Description

11-nor-9-carboxy Cannabinol (Item No. 34592) is an analytical reference standard categorized as a phytocannabinoid metabolite.¹ 11-nor-9-carboxy Cannabinol is a metabolite of cannabinol (Item Nos. ISO60183 | 25495). This product is intended for research and forensic applications.

Reference

1. Watanabe, K., Yamaori, S., Funahashi, T., *et al.* Cytochrome P450 enzymes involved in the metabolism of tetrahydrocannabinols and cannabinol by human hepatic microsomes. *Life Sci.* **80(15)**, 1415-1419 (2007).

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, 01/05/2026

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