

PRODUCT INFORMATION

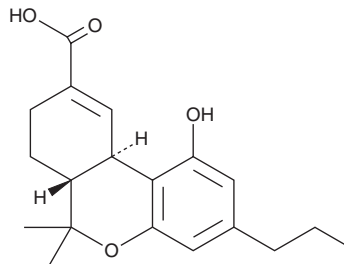


(-)-11-nor-9-carboxy- Δ^9 -Tetrahydrocannabivarin

Item No. 44214

CAS Registry No.: 96912-65-1
Formal Name: 6aR,7,8,10aR-tetrahydro-1-hydroxy-6,6-dimethyl-3-propyl-6H-dibenzo[b,d]pyran-9-carboxylic acid
Synonyms: (-)-11-nor- Δ^9 -Tetrahydrocannabivarin-9-Carboxylic Acid, (-)-11-COOH- Δ^9 -THCV, (-)-11-nor-9-carboxy- Δ^9 -THCV, (-)-11-nor- Δ^9 -THCV-9-COOH

MF: C₁₉H₂₄O₄
FW: 316.4
Purity: ≥98%
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

11-nor-9-carboxy- Δ^9 -Tetrahydrocannabivarin (Item No. 44214) is an analytical reference standard that is structurally similar to known phytocannabinoids. 11-nor-9-carboxy- Δ^9 -Tetrahydrocannabivarin is a potential metabolite of tetrahydrocannabivarin (Item Nos. 18091 | 41381 | 30748) based on the published metabolism of Δ^9 -THC (Item No. 43003 | ISO60157 | 42138).¹ At the time 11-nor-9-carboxy- Δ^9 -tetrahydrocannabivarin (Item No. 44214) was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

Reference

1. Coles, R., Clements, T.T., Nelson, G.J., *et al.* Simultaneous analysis of the Δ^9 -THC metabolites 11-nor-9-carboxy- Δ^9 -THC and 11-hydroxy- Δ^9 -THC in meconium by GC-MS. *J. Anal. Toxicol.* **29(6)**, 522-527 (2016).

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, 10/22/2025

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