

PRODUCT INFORMATION



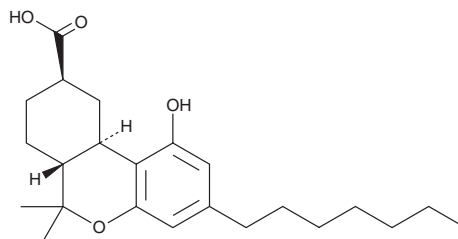
11-nor-9(R)-carboxy-Hexahydrocannabiphorol

Item No. 43122

Formal Name: (6aR,9R,10aR)-3-heptyl-1-hydroxy-6,6-dimethyl-6a,7,8,9,10,10a-hexahydro-6H-benzo[c]chromene-9-carboxylic acid

Synonyms: 11-carboxy-9 β -Hexahydrocannabiphorol, 11-carboxy-HHCP, 11-carboxy-9 β -HHCP, 11-carboxy-9(R)-HHCP, 9(R)-HHCP-COOH, 11-nor-9 β -carboxy-Hexahydrocannabiphorol, 11-nor-9 β -carboxy-HHCP

MF: C₂₃H₃₄O₄
FW: 374.5
Purity: $\geq$ 98%
Supplied as: A neat solid
Storage: -20°C
Stability: $\geq$ 4 years



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

Description

11-nor-9(R)-carboxy-Hexahydrocannabiphorol (Item No. 43122) is an analytical reference standard categorized as a phytocannabinoid metabolite.¹ 11-nor-9(R)-carboxy-Hexahydrocannabiphorol is a metabolite of 9(R)-hexahydrocannabiphorol (Item No. 36346). This product is intended for research and forensic applications.

Reference

1. Lindbom, K., Norman, C., Baginski, S., *et al.* Human metabolism of the semi-synthetic cannabinoids hexahydrocannabinol, hexahydrocannabiphorol and their acetates using hepatocytes and urine samples. *Drug Test Anal.* (2024).

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, 11/26/2025

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