

PRODUCT INFORMATION



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

Item No. 36355

CAS Registry No.: 64663-38-3

Formal Name: [6aR-(6a α ,9 α ,10a β)]-6a,7,8,9,10,10a-hexahydro-1-hydroxy-6,6-dimethyl-3-pentyl-6H-dibenzo[b,d]pyran-9-carboxylic acid

Synonyms: 11-carboxy-9 β -Hexahydrocannabinol, 11-carboxy-HHC, 11-carboxy-9 β -HHC, 11-carboxy-9(R)-HHC, 9(R)-HHC-COOH, 11-nor-9 β -carboxy-Hexahydrocannabinol, 11-nor-9 β -carboxy-HHC

MF: C₂₁H₃₀O₄

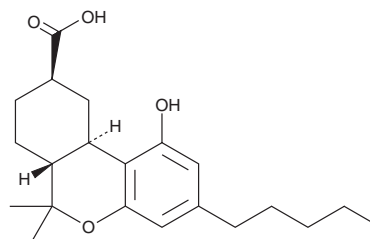
FW: 346.5

Purity: $\geq$ 98%

Supplied as: A solid

Storage: -20°C

Stability: $\geq$ 5 years



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

Description

11-nor-9(R)-carboxy-Hexahydrocannabinol (Item No. 36355) is an analytical reference standard that is structurally similar to known phytocannabinoids.¹ 11-nor-9(R)-carboxy-Hexahydrocannabinol is a metabolite of exo-THC (Item No. 36898) and 9-hexahydrocannabinol.¹⁻³ This product is intended for research and forensic applications.

References

1. Harvey, D.J. The mass spectra of the trimethylsilyl derivatives of *cis*- and *trans*- hexahydrocannabinol and their hydroxy and acid analogues. *Biomed. Mass Spectrom.* **8**(8), 366-372 (1981).
2. Di Trana, A., Di Giorgi, A., Sprega, G., *et al.* Disposition of hexahydrocannabinol epimers and their metabolites in biological matrices following a single administration of smoked hexahydrocannabinol: A preliminary study. *Pharmaceuticals (Basel)* **17**(2), 249 (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, 11/24/2025

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