

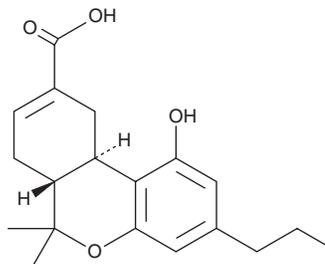
PRODUCT INFORMATION



(-)-11-nor-9-carboxy- Δ^8 -Tetrahydrocannabivarin

Item No. 36631

Formal Name:	1-hydroxy-6,6-dimethyl-3-propyl-6aR,7,10,10aR-tetrahydro-6H-benzo[c]chromene-9-carboxylic acid
Synonym:	(-)-11-nor-9-carboxy- Δ^8 -THCV, (-)-11-nor-9-carboxy- Δ^8 -THV
MF:	C ₁₉ H ₂₄ O ₄
FW:	316.4
Purity:	≥98%
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

(-)-11-nor-9-carboxy- Δ^8 -Tetrahydrocannabivarin (Item No. 36631) is an analytical reference standard that is structurally similar to known phytocannabinoids. (-)-11-nor-9-carboxy- Δ^8 -Tetrahydrocannabivarin is a potential metabolite of Δ^8 -tetrahydrocannabivarin (Item No. 27322) based on the published metabolism of Δ^8 -THC (Item Nos. ISO60158 | 14042).¹ At the time (-)-11-nor-9-carboxy- Δ^8 -tetrahydrocannabivarin (Item No. 36631) 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. Watanabe, K., Yamamoto, I., Oguri, K., *et al.* Comparison in mice of pharmacological effects of Δ^8 -tetrahydrocannabinol and its metabolites oxidized at 11-position. *Eur. J. Pharm.* **63(1)**, 1-6 (1980).

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