

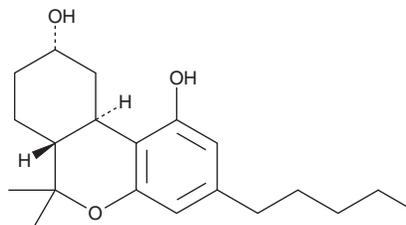
# PRODUCT INFORMATION



## (±)-9-nor-9α-hydroxy Hexahydrocannabinol

Item No. 36256

**CAS Registry No.:** 56469-26-2  
**Formal Name:** *rel*-6aR,7,8,9S,10,10aR-hexahydro-6,6-dimethyl-3-pentyl-6H-dibenzo[b,d]pyran-1,9-diol  
**Synonym:** 9-nor-9α-hydroxy HHC  
**MF:** C<sub>20</sub>H<sub>30</sub>O<sub>3</sub>  
**FW:** 318.5  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

(±)-9-nor-9α-hydroxy Hexahydrocannabinol (Item No. 36256) is an analytical reference standard that is structurally similar to known synthetic cannabinoids.<sup>1</sup> This product is intended for research and forensic applications.

### Reference

1. Wilson, R.S. and May, E.L. 9-Nor-9-hydroxyhexahydrocannabinols. Synthesis, some behavioral and analgesic properties, and comparison with the tetrahydrocannabinols. *J. Med. Chem.* **19(9)**, 1165-1167 (1976).

#### 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, 07/18/2023

#### 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](http://WWW.CAYMANCHEM.COM)