

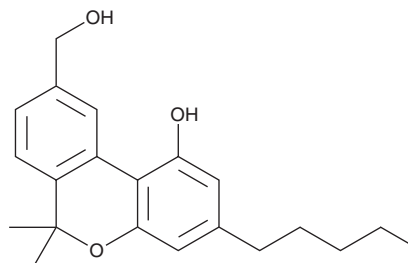
# PRODUCT INFORMATION



## 11-hydroxy Cannabinol

Item No. 34591

**CAS Registry No.:** 30432-08-7  
**Formal Name:** 1-hydroxy-6,6-dimethyl-3-pentyl-6H-dibenzo[b,d]pyran-9-methanol  
**Synonym:** 11-hydroxy CBN  
**MF:** C<sub>21</sub>H<sub>26</sub>O<sub>3</sub>  
**FW:** 326.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224, 286 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Synthetic



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

### Description

11-hydroxy Cannabinol (Item No. 34591) is an analytical reference standard categorized as a phytocannabinoid metabolite.<sup>1-3</sup> 11-hydroxy Cannabinol is an active metabolite of cannabinol (Item Nos. ISO60183 | 25495). This product is intended for research and forensic applications.

### References

1. Yamamoto, I., Watanabe, K., Kuzuoka, K., *et al.* The pharmacological activity of cannabinol and its major metabolite, 11-hydroxycannabinol. *Chem. Pharm. Bull. (Tokyo)* **35(5)**, 2144-2147 (1987).
2. Rhee, M.H., Vogel, Z., Barg, J., *et al.* Cannabinol derivatives: Binding to cannabinoid receptors and inhibition of adenylylcyclase. *J. Med. Chem.* **40(20)**, 3228-3233 (1997).
3. Watanabe, K., Yamaori, S., Funahashi, T., *et al.* Cytochrome P450 enzymes involved in the metabolism of tetrahydrocannabinols and cannabinol by human hepatic microsomes. *Life Sci.* **80(15)**, 1415-1419 (2007).

#### 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, 07/12/2023

#### 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