

# PRODUCT INFORMATION



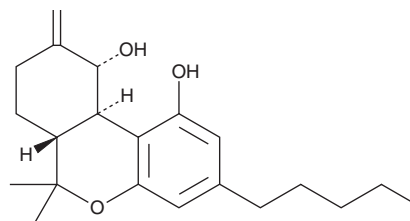
## 10 $\alpha$ -hydroxy-exo-THC

Item No. 38020

CAS Registry No.: 56986-42-6

Formal Name: [6aR-(6 $\alpha$ ,10 $\beta$ ,10 $\alpha$ \beta)]-6a,7,8,9,10,10a-hexahydro-6,6-dimethyl-9-methylene-3-pentyl-6H-dibenzo[b,d]pyran-1,10-diol

Synonyms: 10(R)-hydroxy- $\Delta^9(11)$ -Hexahydrocannabinol, 10 $\alpha$ -hydroxy- $\Delta^9(11)$ -Hexahydrocannabinol, 10 $\alpha$ -hydroxy- $\Delta^9(11)$ -HHC, 10(R)-hydroxy-exo-Tetrahydrocannabinol, 10(R)-hydroxy-exo-THC, 10 $\alpha$ -hydroxy-exo-Tetrahydrocannabinol, 2(R)-hydroxy- $\Delta^1(7)$ -Tetrahydrocannabinol, 2 $\alpha$ -hydroxy- $\Delta^1(7)$ -Tetrahydrocannabinol, 10 $\alpha$ -hydroxy- $\Delta^9(11)$ -Tetrahydrocannabinol, 2(R)-hydroxy- $\Delta^1(7)$ -THC, 2 $\alpha$ -hydroxy- $\Delta^1(7)$ -THC, 10 $\alpha$ -hydroxy- $\Delta^9(11)$ -THC



MF: C<sub>21</sub>H<sub>30</sub>O<sub>3</sub>

FW: 330.5

Purity:  $\geq 98\%$

UV/Vis.:  $\lambda_{\max}$ : 210 nm

Supplied as: A 10 mg/ml solution in acetonitrile

Storage: -20°C

Stability:  $\geq 2$  years

Item Origin: Synthetic

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

### Description

10 $\alpha$ -hydroxy-exo-THC (Item No. 38020) is an analytical reference standard categorized as a phytocannabinoid.<sup>1</sup> 10 $\alpha$ -hydroxy-exo-THC is an active metabolite of  $\Delta^9$ -THC (Item Nos. ISO60157 | 12068).<sup>1,2</sup> It has been found in *Cannabis*.<sup>1</sup> 10 $\alpha$ -hydroxy-exo-THC is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

### References

1. Radwan, M.M., ElSohly, M.A., El-Alfry, A.T., *et al.* Isolation and pharmacological evaluation of minor cannabinoids from high-potency *Cannabis sativa*. *J. Nat. Prod.* **78(6)**, 1271-1276 (2015).
2. Pitt, C.G., Fowler, M.S., Sathe, S., *et al.* Synthesis of metabolites of  $\Delta^9$ -tetrahydrocannabinol. *J. Am. Chem. Soc.* **97(13)**, 3798-3802 (1975).

#### 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, 11/11/2025

#### 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