

PRODUCT INFORMATION



10(S)-hydroxy-9(R)-Hexahydrocannabinol

Item No. 38273

CAS Registry No.: 73648-83-6

Formal Name: 6aR-6a α ,7,8,9 α ,10 β ,10a β -hexahydro-6,6,9-trimethyl-3-pentyl-6H-dibenzo[b,d]pyran-1,10-diol

Synonyms: 10 α -hydroxy-9(R)-Hexahydrocannabinol, 10 α -hydroxy-9(R)-HHC, 10(S)-hydroxy-9(R)-HHC

MF: C₂₁H₃₂O₃

FW: 332.5

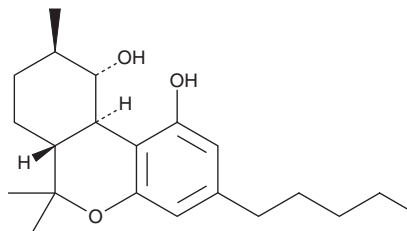
Purity: $\geq 98\%$

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

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

10(S)-hydroxy-9(R)-Hexahydrocannabinol (Item No. 38273) is an analytical reference standard that is structurally similar to known phytocannabinoids. 10(S)-hydroxy-9(R)-Hexahydrocannabinol has been found in Cannabis.¹ It induces drowsiness, hypolocomotion, and occasional partial ptosis and head drop, markers of psychotropic activity, in rhesus monkeys.² 10(S)-hydroxy-9(R)-Hexahydrocannabinol is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Ahmed, S.A., Ross, S.A., Slade, D., *et al.* Minor oxygenated cannabinoids from high potency *Cannabis sativa* L. *Phytochemistry* **117**, 194-199 (2015).
2. Mechoulam, R., Lander, N., Varkony, T.H., *et al.* Stereochemical requirements for cannabinoid activity. *J. Med. Chem.* **23**(10), 1068-1072 (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, 01/06/2026

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