

# Hexahydrocannabinols (HHCs)

Hydrogenation of tetrahydrocannabinols ( $\Delta^8$ - or  $\Delta^9$ -THC) leads to formation of a mixture of hexahydrocannabinols (HHCs) comprised of the 9(S)- and 9(R)-HHC diastereomers. These compounds retain some psychoactivity but avoid classification as THC as well as any THC-related regulations. Separation of the HHC diastereomers is possible by chromatography (HPLC, GC), but identity can only be confirmed with verified reference standards. Cayman offers fully characterized reference standards to aid in the correct identification and differentiation of hydrogenated phytocannabinoids.



ACCREDITED  
ISO/IEC 17025 #AT-1773  
ISO 17034 #AR-1774

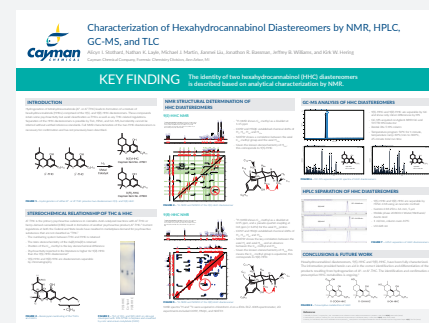
## Explore Cayman's Scientific Poster

### Characterization of Hexahydrocannabinol Diastereomers by NMR, HPLC, GC-MS, and TLC

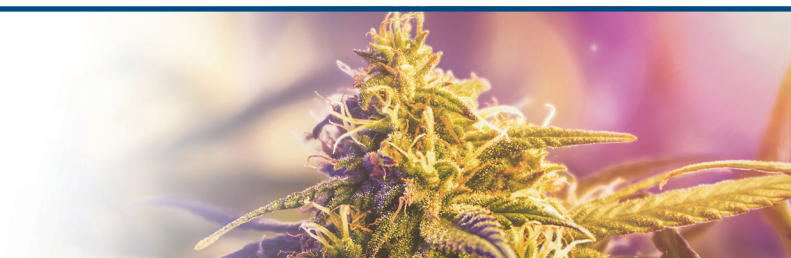
Alicyn I. Stothard, Nathan K. Layle, Michael J. Martin, Jianmei Liu, Jonathon R. Bassman, Jeffrey B. Williams, and Kirk W. Hering - Cayman Chemical Forensic Division

Discover a reliable method to differentiate two HHC diastereomers resulting from  $\Delta^8$ - or  $\Delta^9$ -THC hydrogenation.

[View the full poster inside](#)



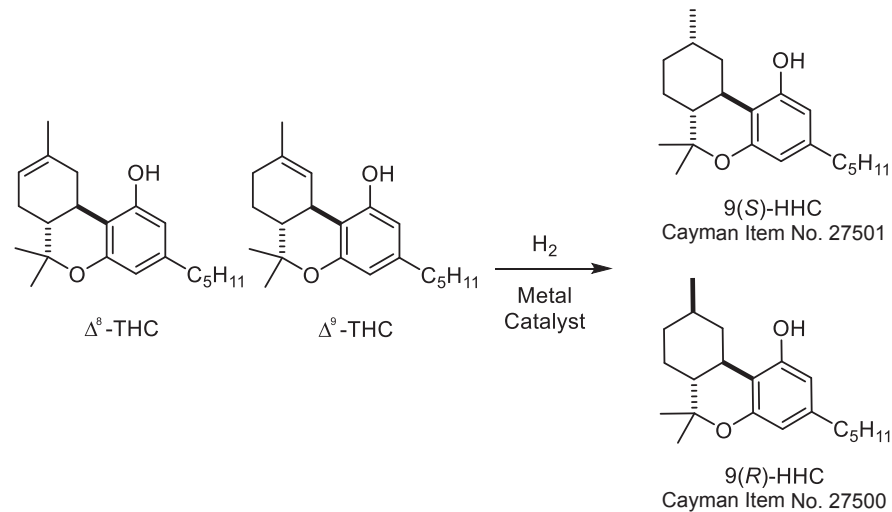
Discover all  
HHC Cannabinoid  
Analytical Standards



**KEY FINDING** The identity of two hexahydrocannabinol (HHC) diastereomers is described based on analytical characterization by NMR.

## INTRODUCTION

Hydrogenation of tetrahydrocannabinols ( $\Delta^8$ - or  $\Delta^9$ -THC) leads to formation of a mixture of hexahydrocannabinols (HHCs) comprised of the 9(S)- and 9(R)-HHC diastereomers. These compounds retain some psychoactivity but avoid classification as THC<sub>s</sub> as well as any THC-related regulations. Separation of the HHC diastereomers is possible by TLC, HPLC, and GC-MS, but identity cannot be inferred without verified reference standards. Full NMR characterization of the two HHC diastereomers is necessary for confirmation and has not previously been described.

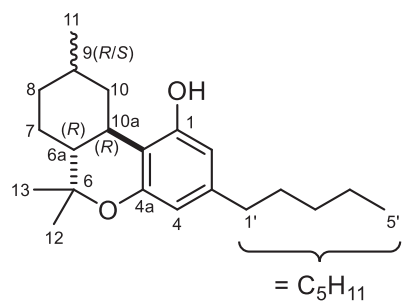


**FIGURE 1** – Hydrogenation of either  $\Delta^8$ - or  $\Delta^9$ -THC provides two diastereomers 9(S)- and 9(R)-HHC

## STEREOCHEMICAL RELATIONSHIP OF THC & HHC

$\Delta^9$ -THC is the primary psychoactive substance in *Cannabis*. Acid-catalyzed reactions with  $\Delta^9$ -THC or hemp-derived cannabidiol (CBD) result in formation of another psychoactive product,  $\Delta^8$ -THC.<sup>1</sup> Current regulations at both the Federal and State levels have resulted in marketplace demand for psychoactive substances that are not classified as “THC.”

- The numbering system between THC and HHC is retained
- The *trans* stereochemistry of the 6a(R),10a(R) is retained
- Position of the C<sub>11</sub>-methyl is the key stereochemical difference
- Psychoactivity reported in the literature is higher for 9(R)-HHC than the 9(S)-HHC diastereomer<sup>2</sup>
- 9(S)-HHC and 9(R)-HHC are diastereomers separable by chromatography



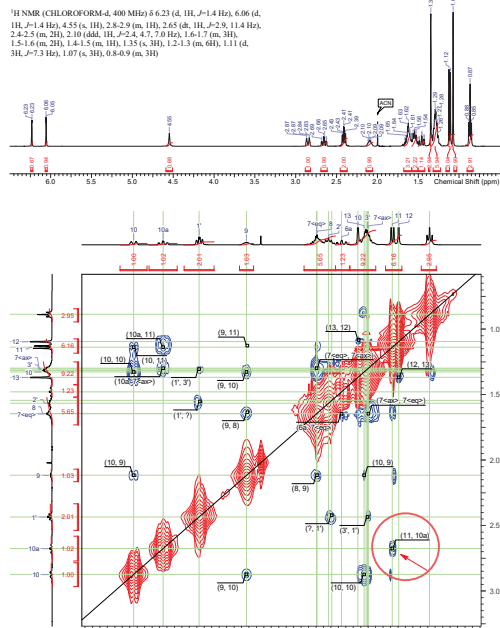
**FIGURE 2** – Benzopyran numbering of the THC<sub>s</sub> and HHC<sub>s</sub>



**FIGURE 3** – TLC of 9(S)- and 9(R)-HHC on silica gel developed with 10% MTBE in heptane and visualized by ceric ammonium molybdate (CAM)

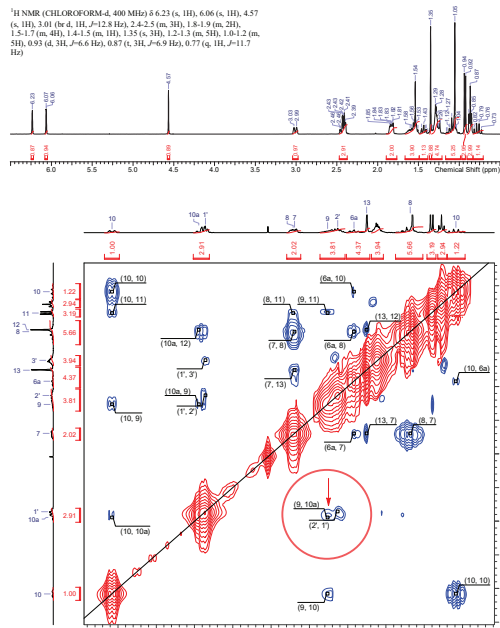
## NMR STRUCTURAL DETERMINATION OF HHC DIASTEREOMERS

## 9(S)-HHC NMR



**FIGURE 4** –  $^1\text{H}$ -NMR and NOESY of the 9(*S*)-HHC diastereomer

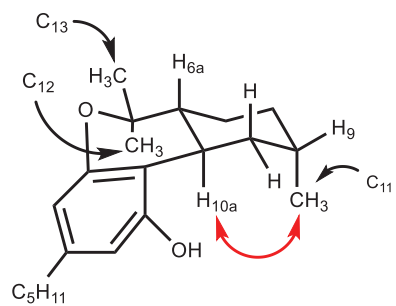
## 9(R)-HHC NMR



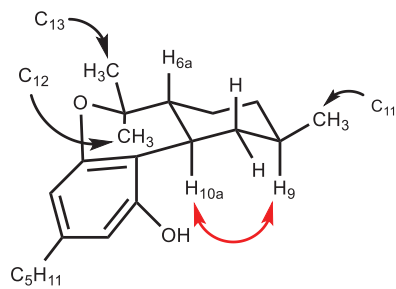
**FIGURE 5** –  $^1\text{H}$ -NMR and NOESY of the 9(*R*)-HHC diastereomer

NMR spectra  $^1\text{H}$  and  $^{13}\text{C}$  were acquired in chloroform-d on a JEOL ECZ-400S spectrometer, 2D experiments included COSY, HSQC, and NOESY.

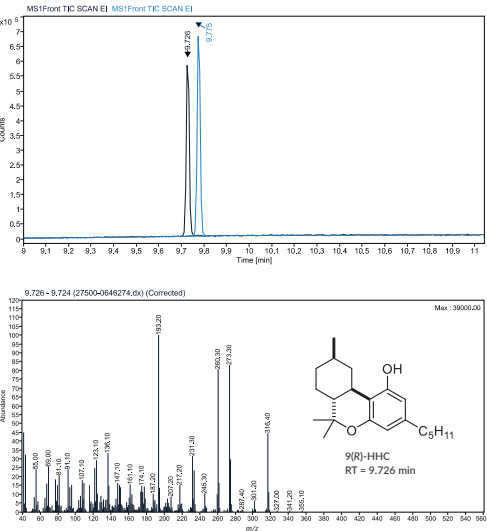
- $^1\text{H-NMR}$  shows  $\text{C}_{11}$ -methyl as a doublet at 1.15 ppm
- COSY and HSQC established chemical shifts of  $\text{H}_{9'}$ ,  $\text{H}_{10'}$ ,  $\text{H}_{6a'}$  and  $\text{H}_{10a}$
- NOESY shows a correlation between the axial  $\text{C}_{11}$ -methyl group and the axial  $\text{H}_{10a}$
- Given the known stereochemistry of  $\text{H}_{10a}$ , this corresponds to 9(*S*)-HHC



- $^1\text{H-NMR}$  shows  $\text{C}_{11}$ -methyl as a doublet at 0.95 ppm, and a pseudo quartet coupling at 0.8 ppm ( $J=12\text{Hz}$ ) for the axial  $\text{H}_{10}$  proton
- COSY and HSQC established chemical shifts of  $\text{H}_9$ ,  $\text{H}_{10}$ ,  $\text{H}_{6a}$  and  $\text{H}_{10a}$
- NOESY shows the key correlation between the axial  $\text{H}_9$  and axial  $\text{H}_{10a}$ , and an absence between the  $\text{C}_{11}$ -methyl and  $\text{H}_{10a}$
- Given the known stereochemistry of  $\text{H}_{10a}$ , this means the  $\text{C}_{11}$ -methyl group is equatorial, this corresponds to 9(R)-HHC



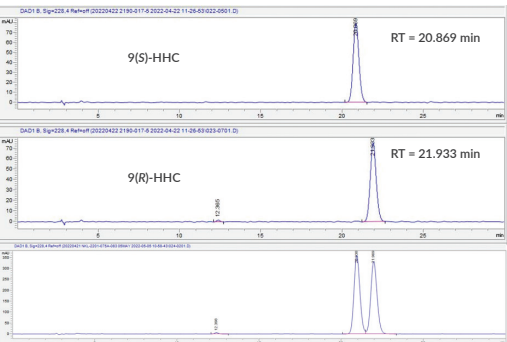
## GC-MS ANALYSIS OF HHC DIASTEREOMERS



**FIGURE 6** – GC-MS separation and EI spectra of HHC diastereomers

- 9(S)-HHC and 9(R)-HHC are separable by GC and show only minor differences by MS
- GC-MS acquired on Agilent 8890 GC and 5977B MS Detector
- Restek Rtx-5 MS column
- Temperature program 50°C for 1 minute, temperature ramp 30°C/min to 300°C, 25-minute total run time

## HPLC SEPARATION OF HHC DIASTEREOMERS

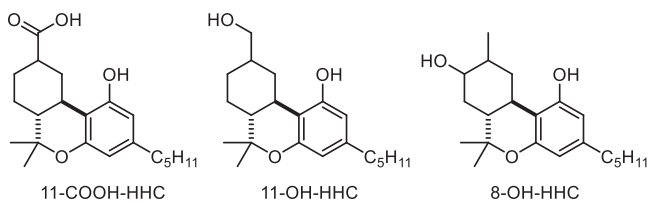


**FIGURE 7** – HPLC separation of HHC diastereomers

- 9(S)-HHC and 9(R)-HHC are separable by HPLC-C18 using an isocratic method
- Gemini-C18 250 x 4.6 mm, 5  $\mu$ m
- Mobile phase 20:80:0.1 Water/Methanol/Acetic Acid
- 1 ml/min, column oven 40°C
- UV 228 nm

## CONCLUSIONS & FUTURE WORK

Hexahydrocannabinol diastereomers, 9(*S*)-HHC and 9(*R*)-HHC, have been fully characterized. The information provided herein can aid in the correct identification and differentiation of the products resulting from hydrogenation of  $\Delta^8$ - or  $\Delta^9$ -THC. The identification and confirmation of presumptive HHC metabolites is ongoing.<sup>3</sup>



**FIGURE 8 – Presumptive metabolites of HHC**

## References

1. Marzullo, P., Foschi, F., Coppini, D.A., et al. Cannabidiol as the substrate in acid-catalyzed intramolecular cyclization. *J. Nat. Prod.* **83**(10), 2894–2901 (2020).
2. Mechoulam, R., Lander, N., Varkony, T.H., et al. Stereochemical requirements for cannabinoid activity. *J. Med. Chem.* **23**(10), 1068–1072 (1980).
3. Harvey, D.J. and Brown, N.K. In vitro metabolism of the equatorial C<sub>1</sub>-methyl isomer of hexahydrocannabinol in several mammalian species. *Drug Metab. Dispos.* **19**(3), 714–716 (1991).

# Hydrogenated Phytocannabinoids Available from Cayman

More than 20 reference standards for:

- Hexahydrocannabinols (HHCs)
- Hexahydrocannabiphorols (HHCPs)
- Tetrahydrocannabinidiols (H4-CBDs)

## Hexahydrocannabinols (HHCs)

Item No.	Product Name
27500	9(R)-Hexahydrocannabinol
37913	9(R)-Hexahydrocannabinol (CRM)
35368	9(R)-Hexahydrocannabinol Acetate
27501	9(S)-Hexahydrocannabinol
37942	9(S)-Hexahydrocannabinol (CRM)
35369	9(S)-Hexahydrocannabinol Acetate

## Other HHCs of Interest

Item No.	Product Name
36256	(±)-9-nor-9 $\alpha$ -hydroxy Hexahydrocannabinol
36257	(±)-9-nor-9 $\beta$ -hydroxy Hexahydrocannabinol
36129	(±)-9 $\alpha$ -hydroxy Hexahydrocannabinol
35266	(±)-9 $\beta$ -hydroxy Hexahydrocannabinol
36353	11-hydroxy-9(R)-Hexahydrocannabinol
36354	11-hydroxy-9(S)-Hexahydrocannabinol
36355	11-nor-9(R)-carboxy-Hexahydrocannabinol
36356	11-nor-9(S)-carboxy-Hexahydrocannabinol
36250	8(R)-hydroxy-9(R)-Hexahydrocannabinol
36249	8(S)-hydroxy-9(S)-Hexahydrocannabinol

View all HHC cannabinoids at [www.caymanchem.com](http://www.caymanchem.com)

## Hexahydrocannabiphorols (HHCPs)

Item No.	Product Name
36346	9(R)-Hexahydrocannabiphorol
37847	9(R)-Hexahydrocannabiphorol Acetate
36347	9(S)-Hexahydrocannabiphorol
37848	9(S)-Hexahydrocannabiphorol Acetate

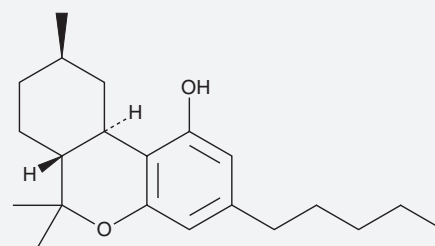
View all HHCP cannabinoids at [www.caymanchem.com](http://www.caymanchem.com)

## Tetrahydrocannabinidiols (H4-CBDs)

Item No.	Product Name
36350	1(R)-Tetrahydrocannabinidiol (1(R)-H4-CBD)
36351	1(S)-Tetrahydrocannabinidiol (1(S)-H4-CBD)

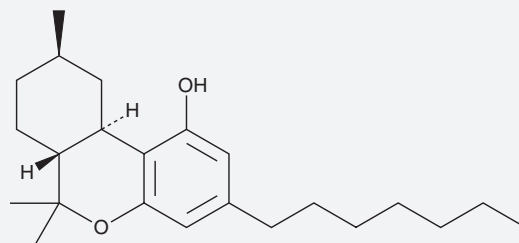
View all H4-CBD cannabinoids at [www.caymanchem.com](http://www.caymanchem.com)

## Major Hydrogenated Phytocannabinoids



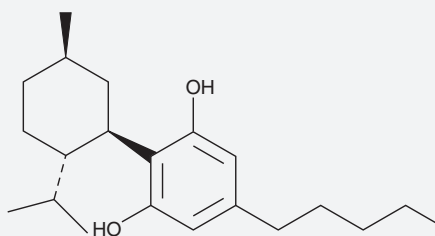
**9(R)-HHC**

Item No. 27500



**9(R)-HHCP**

Item No. 36346



**1(R)-H4-CBD**

Item No. 36350

## Can't find your unknown?

Submit your GC-MS data (preferred) and any relevant information (suspected chemical class, matrix, solvent, GC parameters, etc.) to [techserv@caymanchem.com](mailto:techserv@caymanchem.com) and our scientists can help you solve your unknown.

**Visit our Cannabis & Hemp  
Resource Center  
to learn more**