

## Application Note

# GC Separation for Identification of *iso*-THC Contaminants and Accurate Quantification of $\Delta^8$ -THC and $\Delta^9$ -THC in *Cannabis* Samples

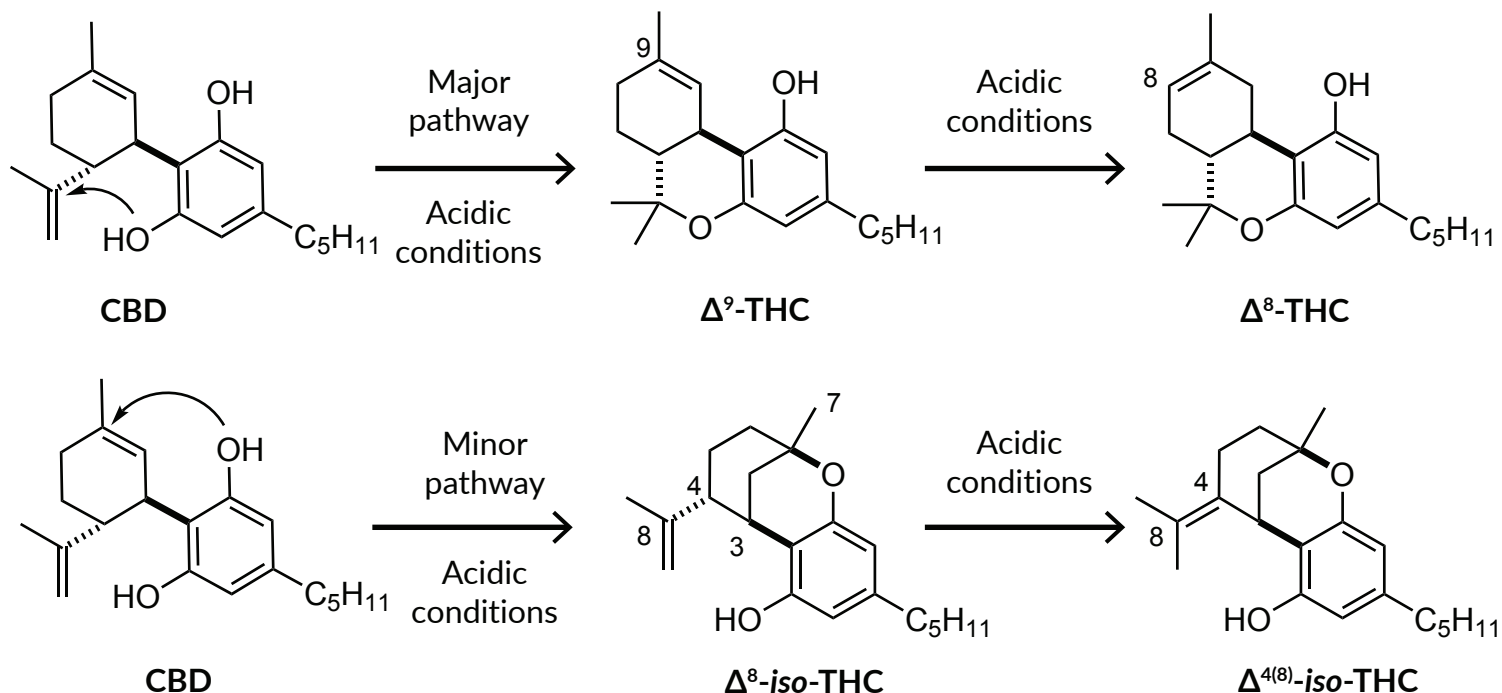
Jeffrey B. Williams, Jianmei Liu, and Kirk W. Hering, Ph.D.  
Cayman Chemical Company, Ann Arbor, MI

## Key Features

- Synthetic conversion of cannabidiol (CBD) to the tetrahydrocannabinols (THCs)  $\Delta^9$ -THC and  $\Delta^8$ -THC produces measurable quantities of the *iso*-THC products  $\Delta^8$ -*iso*-THC and  $\Delta^{4(8)}$ -*iso*-THC.
- Co-elution of THCs and *iso*-THCs makes C18 reversed phase-HPLC (RP-HPLC) methods unsuitable for the quantification of THCs. Quantities of THCs are distorted if the testing methods are unable to resolve the *iso*-THCs in analyzed samples.
- Gas chromatography (GC) offers a reliable and robust method for quantification of  $\Delta^9$ -THC,  $\Delta^8$ -THC,  $\Delta^8$ -*iso*-THC, and  $\Delta^{4(8)}$ -*iso*-THC. Additionally, this method represents a method for identifying THC derived from CBD conversion.

# Introduction

Analytical testing of *Cannabis* products for identification and accurate quantification of psychoactive components such as  $\Delta^9$ -THC and  $\Delta^8$ -THC is a regulatory requirement for safety and distribution. However, quantification of THC<sub>s</sub> may be complicated by the introduction of the closely eluting side products  $\Delta^8$ -*iso*-THC and  $\Delta^{4(8)}$ -*iso*-THC.<sup>1-3</sup> These constitutional isomers of THC form during synthetic conversion of CBD under acidic conditions (Figure 1).



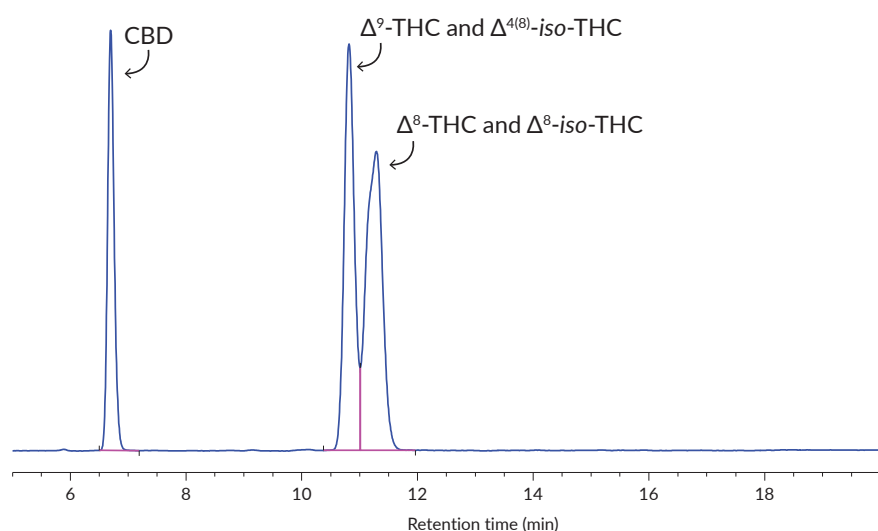
**Figure 1.** Acid-catalyzed cyclization of CBD to form THC<sub>s</sub> and *iso*-THC<sub>s</sub>.

The *iso*-THC<sub>s</sub> were first identified by Mechoulam's research group in a 1968 paper where CBD treated with the Lewis acid boron-trifluoride diethyl etherate led to formation of THC with 13% of  $\Delta^8$ -*iso*-THC isolated as a side product. Mechoulam also identified  $\Delta^{4(8)}$ -*iso*-THC in this same paper and demonstrated that it could be formed by boiling  $\Delta^8$ -*iso*-THC in a mixture of sulfuric acid and methanol.<sup>4</sup> A 2020 *Journal of Natural Products* publication by the Passarella group also highlights the formation of these two isomers under a number of acidic conditions and provides thorough NMR characterization.<sup>5</sup> To date, the *iso*-THC<sub>s</sub> have not been tested for pharmacological activity or safety in any published results. Since the *iso*-THC<sub>s</sub> have not been found naturally in any *Cannabis sativa* L., their presence in *Cannabis* products serves as a useful marker to identify THC derived from the synthetic conversion of CBD.

This application note highlights the co-elution observed between THC<sub>s</sub> and *iso*-THC<sub>s</sub> by RP-HPLC and presents a GC method that provides for baseline separation and quantification of these species.

## RP-HPLC Methods

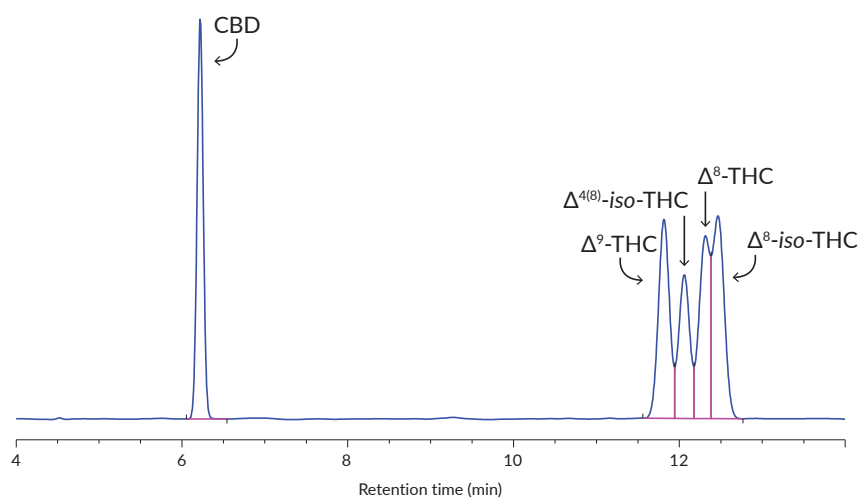
A mixture of equal parts CBD,  $\Delta^9$ -THC,  $\Delta^8$ -THC,  $\Delta^8$ -*iso*-THC, and  $\Delta^{4(8)}$ -*iso*-THC was injected on a Gemini® C18 column using isocratic RP conditions, resulting in significant co-elution of the THC<sub>s</sub> and *iso*-THC<sub>s</sub> (**Figure 2**). In conventional samples of CBD-derived THC<sub>s</sub>, the amount of *iso*-THC<sub>s</sub> present is likely to be a much smaller percentage relative to the THC<sub>s</sub> and would almost certainly be hidden underneath the THC peaks with similar chromatographic conditions. Consequently, the presence of these side products would likely be presented as a false enhancement of the co-eluting THC peak.



Analyte	Retention time (min)	Item No.
CBD	6.62	ISO60156
$\Delta^9$ -THC	10.828	ISO60157
$\Delta^{4(8)}$ - <i>iso</i> -THC	10.828	33863
$\Delta^8$ -THC	11.115	ISO60158
$\Delta^8$ - <i>iso</i> -THC	11.287	33864

**Figure 2.** Co-injection of CBD,  $\Delta^9$ -THC,  $\Delta^{4(8)}$ -*iso*-THC,  $\Delta^8$ -THC, and  $\Delta^8$ -*iso*-THC on a Gemini® C18 column under RP conditions. Data was acquired on an Agilent 1100 Series HPLC with a Gemini® C18 column (250 mm x 4.6 mm, 5  $\mu$ m) using a 20:80 water:acetonitrile with 0.1% acetic acid mobile phase, 1 ml/min flow rate, 40°C column temperature, and monitoring UV at 228 nm.

Efforts to find a more optimized RP-HPLC method involved screening several columns and conditions. Improved results were observed using a Raptor® ARC-C18 HPLC column under isocratic RP conditions (**Figure 3**). However, this method still results in significant overlap of the *iso*-THC and THC analytes and a lack of baseline resolution. As such, GC was explored next as another possible way to fully resolve these isomers.

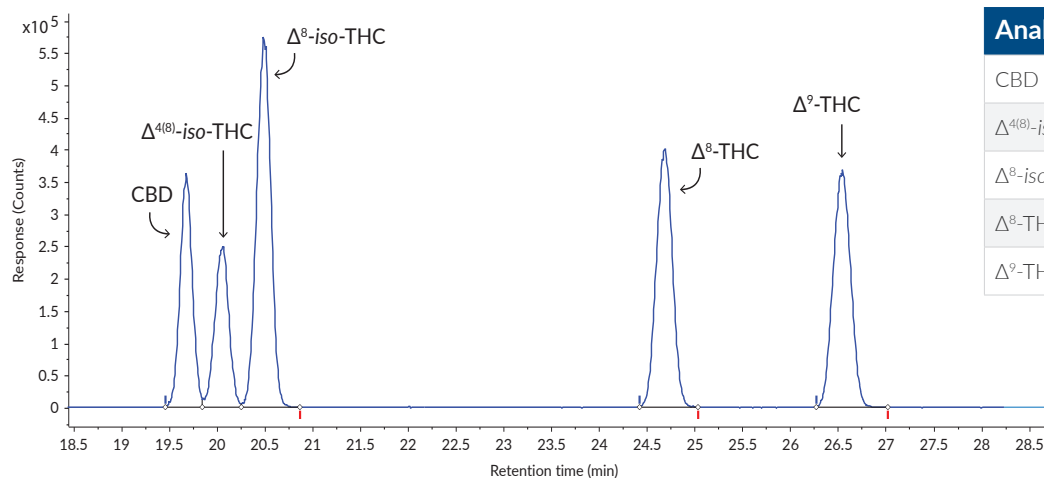


Analyte	Retention time (min)	Item No.
CBD	6.221	ISO60156
$\Delta^9$ -THC	11.816	ISO60157
$\Delta^{4(8)}$ - <i>iso</i> -THC	12.062	33863
$\Delta^8$ -THC	12.317	ISO60158
$\Delta^8$ - <i>iso</i> -THC	12.469	33864

**Figure 3.** Co-injection of CBD,  $\Delta^9$ -THC,  $\Delta^{4(8)}$ -*iso*-THC,  $\Delta^8$ -THC, and  $\Delta^8$ -*iso*-THC on a Raptor® ARC-C18 column under RP conditions. Data was acquired on an Agilent 1100 Series HPLC with a Raptor® ARC-C18 column (150 mm x 4.6 mm, 2.7  $\mu$ m) using a 30:70 water:acetonitrile with 0.1% acetic acid mobile phase, 1 ml/min flow rate, 40°C column temperature, and monitoring UV at 228 nm.

# GC Method

An optimized GC method for separating a mixture of CBD,  $\Delta^9$ -THC,  $\Delta^8$ -THC,  $\Delta^8$ -iso-THC, and  $\Delta^{4(8)}$ -iso-THC resulted in baseline resolution of all the peaks and is shown in **Figure 4**. One of the highlights of this method is that the iso-THCs elute closer to CBD and have greater than four minutes separation in retention time from the THC. Hence, this provides an excellent means to determine the quantity of iso-THCs present in a sample with complete resolution from the THCs. While this method was demonstrated on a GC-MS instrument, the conditions may be amended for GC-FID instruments.



Analyte	Retention time (min)	Item No.
CBD	19.675	ISO60156
$\Delta^{4(8)}$ -iso-THC	20.066	33863
$\Delta^8$ -iso-THC	20.482	33864
$\Delta^8$ -THC	24.683	ISO60158
$\Delta^9$ -THC	26.546	ISO60157

**Figure 4.** Co-injection of CBD,  $\Delta^{4(8)}$ -iso-THC,  $\Delta^8$ -iso-THC,  $\Delta^8$ -THC, and  $\Delta^9$ -THC on GC method.

Data was acquired on an Agilent 8890 GC equipped with a 5977B MS detector, using a Restek® Rtx-5MS capillary column (30 m x 320  $\mu$ m x 0.5  $\mu$ m). The injector temperature was set at 300°C, and the oven temperature was programmed to start at 50°C and increase 40°C/min to 210°C, hold 20 min, then increase 40°C/min to 300°C and hold 8.75 min. The total run time was 35 min. The injection mode was set to split with a ratio of 60:1. Helium was used as the carrier gas at a constant flow of 2 ml/min.

## Conclusion

Quantification of  $\Delta^9$ -THC and  $\Delta^8$ -THC content in extracts, edibles, and other *Cannabis*-derived products has become increasingly important as these materials are legalized for medicinal and recreational use. Both synthetic production and chemical conversion of plant-derived materials may result in the formation of iso-THCs, closely eluting side products of THCs. While most RP-HPLC methods may not fully resolve the iso-THCs from  $\Delta^9$ -THC and  $\Delta^8$ -THC, we present here a GC method that can be used to fully resolve these compounds, aiding in the development of testing methods to identify and accurately quantify THCs and iso-THCs in *Cannabis* products.

## Acknowledgements

Special thanks to Dr. Richard Sams at KCA Laboratories for suggesting the GC method.

## References

1. Meehan-Atrash, J. and Rahman, I. Novel  $\Delta^8$ -tetrahydrocannabinol vaporizers contain unlabeled adulterants, unintended byproducts of chemical synthesis, and heavy metals. *Chem. Res. Toxicol.* **35**(1), 73-76 (2022).
2. Radwan, M.M., Wanas, A.S., Gul, W., et. al. Isolation of impurities in commercially marketed  $\Delta^8$ -THC products. *J. Nat. Prod.* (2023).
3. Geci, M., Scialdone, M., and Tishler, J. The dark side of cannabidiol: The unanticipated social and clinical implications of synthetic  $\Delta^8$ -THC. *Cannabis Cannabinoid Res.* **8**(2), 270-282 (2023).
4. Gaoni, Y. and Mechoulam, R. The iso-Tetrahydrocannabinols. *Israel J. Chem.* **6**(5), 679-690 (1968).
5. 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).