

Fully Automated Sample Preparation for the Analysis of Terpenes in Cannabis Flower

Using the Agilent PAL3 Series II RTC Automation with
an Agilent Intuvo 9000/5977C GC/MS

Authors

Saurabh U. Patel,
Jessica L. Westland,
and Samuel P. Haddad
Agilent Technologies, Inc.

Abstract

Analysis of terpenes in cannabis by gas chromatography/mass spectrometry (GC/MS) has become an important tool for the characterization of taste and smell profiles of commercially available cannabis strains. This application brief demonstrates the capability of the Agilent PAL3 Series II robotic tool change (RTC) sampler to fully automate the sample preparation process for terpenes in cannabis flower. The use of automated sample preparation helps reduce user errors and the amount of sample and solvent needed for analysis. Coupled with an Agilent Intuvo 9000/5977C GC/MS system, PAL3 Series II RTC automation can save time and cut down on energy consumption, leading to a greener high-throughput laboratory.

Introduction

Terpenes are a class of compounds that are responsible for the unique flavor and aroma in each strain of *Cannabis* spp. Analysis of terpenes in cannabis by gas chromatography/mass spectrometry (GC/MS) has become an important tool for the characterization of taste and smell profiles of commercially available cannabis strains. The manual solvent extraction process, while straightforward, can introduce extraction errors at the bench. This application brief demonstrates the capability of the Agilent PAL3 Series II RTC sampler to fully automate the sample preparation process for terpenes in cannabis flower. The PAL3 Series II RTC, paired with the Agilent Intuvo 9000 GC and 5977C MS detector (MSD), creates an easy and robust solution for any cannabis lab looking to analyze terpenes.

Experimental

Chemicals and reagents

Terpenes standard mixes, CAN-TERP-MIX1H and CAN-TERP-MIX2H, each containing 21 terpenes, were purchased from SPEX CertiPrep (Metuchen, NJ, US). The internal standard (ISTD), 2-fluorobiphenyl and ethyl acetate (purity 99.9%) were purchased from Sigma-Aldrich (St. Louis, MO, US). The test matrix for calibration, cold pressed hemp seed oil (blank matrix), was obtained from Nutiva (Richmond, CA, US). Various strains of ground cannabis flower were obtained from the University of Mississippi.

Instrumentation

The GC/MS instrument method parameters followed the methodology described in Agilent application note 5994-2032EN.¹ GC instrument parameters can be found in Table 1, MS instrument parameters in Table 2, and PAL3 Series II RTC instrument configuration in Table 3. Table 4 contains a list of consumable items used for the current application.

Table 1. GC conditions for terpenes analysis.

| Agilent Intuvo 9000 GC, Auto Injector, and Tray | |
|---|---|
| Injection Volume | 1.0 µL |
| Inlet | EPC split/splitless |
| Mode | Split |
| Split Ratio | 150:1 |
| Septum Purge Flow Mode | Standard, 3 mL/min |
| Inlet Temperature | 250 °C |
| Guard Chip Temperature | 250 °C (Track Oven Off) |
| Bus Temperature | 260 °C |
| Oven | Initial: 75 °C (1 min hold) Ramp 1: 5 °C/min to 165 °C Ramp 2: 175 °C/min to 250 °C (10.514 min hold) |
| Column 1 | Agilent J&W DB-Select 624, 30 m × 0.25 mm, 1.4 µm |
| Control Mode, Flow | Constant flow, 2.0 mL/min |
| Column 2 | Agilent J&W DB-Select 624, 30 m × 0.25 mm, 1.4 µm |
| Control Mode, Flow | Constant flow, 2.2 mL/min |
| PSD Purge Flow | 3 mL/min |
| Agilent Intuvo 9000 GC Backflush Parameters | |
| Inlet Pressure (Backflushing) | 2 psi |
| Backflush Pressure | 65 psi |
| Void Volumes | 5 |
| Backflush Time | 3.236 |

Table 2. MS conditions for terpenes analysis.

| Agilent 5977C MSD | |
|---------------------------|-------------------|
| Source | Agilent Extractor |
| Extractor Lens | 9 mm |
| Transfer Line Temperature | 260 °C |
| Source Temperature | 300 °C |
| Quadrupole Temperature | 200 °C |
| Mode | SIM |
| EM Voltage Gain | Variable |
| Solvent Delay | 13 min |
| Tune File | atune.u |

Table 3. PAL 3 configuration for terpenes analysis.

| PAL3 Series II RTC Rail | |
|-------------------------|--|
| Tools | Liquid handling tools |
| Syringes | 10 µL (part number 8010-1307) 25 µL (part number 8010-1310) 1 mL (part number 8010-1326) |
| Control System | PAL Method Composer v1.4 |
| PAL Modules | Vortex mixer, tray holder, large wash module |
| Optional Modules | Fast wash station, dilutor module |

Table 4. Agilent consumables and part numbers used in the method for terpenes analysis.

| Consumable | Description | Part Number |
|-------------------|---|----------------|
| Inlet Septum | Advanced Green septum, nonstick, 11 mm | 5183-4759 |
| Inlet Liner | Universal Ultra Inert low pressure drop inlet liner | 5190-2295 |
| Guard Chip | Agilent Intuvo S/SL Guard Chip | G4587-60565 |
| Column | DB-Select 624 30 m × 250 µm, 1.4 µm (quantity: 2) | 122-0334UI-INT |
| Gaskets | Intuvo polyimide gasket | 5190-9072 |
| Compression Bolts | Intuvo compression bolts | G4581-60260 |
| Detector Tail | Intuvo MS tail | G4590-60009 |
| Extractor Lens | Extractor EI 9 mm lens | G3870-20449 |
| PAL Vials | Screw top, clear, round bottom, 10 mL | 5188-5392 |
| PAL Screw Caps | Steel, magnetic cap, PTFE/silicone septa 18 mm | 5188-2759 |

Automated calibration and sample preparation

All automated preparation steps were scripted using PAL Method Composer (G7388A). Figure 1 details the working instructions for the automated preparation of the calibration curve levels, while Figure 2 details the working instructions for the automated sample preparation.

To evaluate the automated preparation of calibration curves, three separate curves were created using the PAL3 Series II RTC. A total of eight calibration points were used, ranging from 3.83 µg/mL to 490.2 µg/mL. Accuracy and precision were calculated for each terpene (n = 3) using a low (3.83 µg/mL), a mid (30.64 µg/mL), and a high (245.10 µg/mL) calibration level standard.

To evaluate the automated preparation of samples, homogenized cannabis flower was extracted. Two different cannabis flower samples were extracted in quadruplicate (n = 4) to show precision (Figure 3), and six additional cannabis flower samples were extracted (n = 1) to show sample variation (Figure 4).

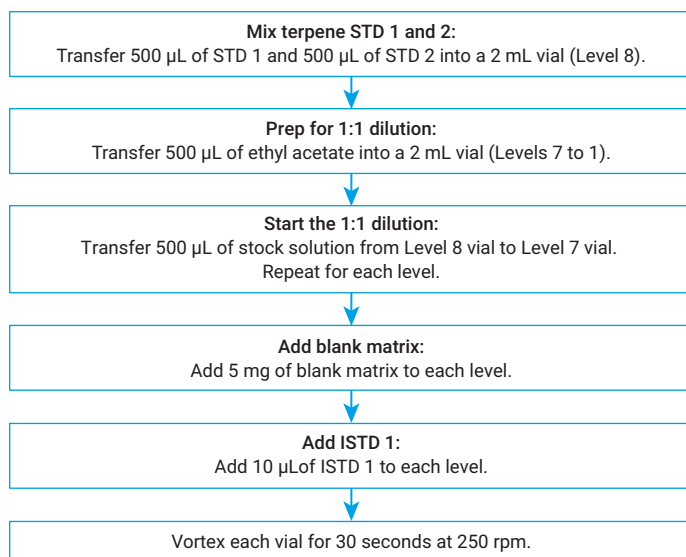


Figure 1. PAL Method Composer working instructions for the automated preparation of terpene calibration curves. The volume needed to accurately transfer 5 mg of hemp seed oil was determined experimentally, while keeping its density and viscosity in mind. Preparation instructions for ISTD 1 (10,000 µg/mL of 2-fluorobiphenyl in ethyl acetate) can be found in Agilent application note 5994-2032EN.¹

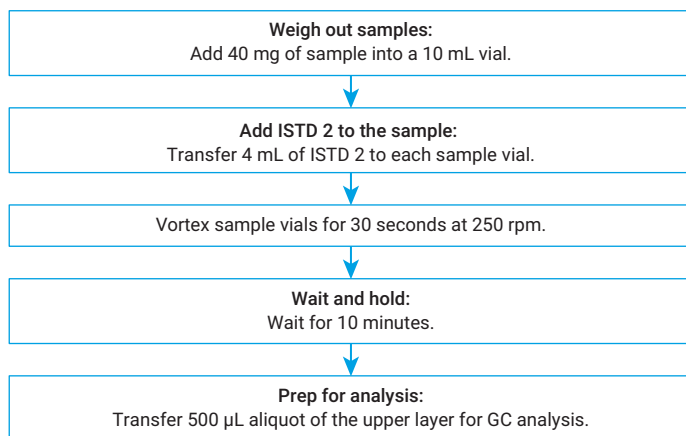


Figure 2. PAL Method Composer working instructions for the automated preparation of terpenes samples. Preparation instruction for ISTD 2 (200 µg/mL of 2-fluorobiphenyl in ethyl acetate) can be found in Agilent application note 5994-2032EN.¹



Figure 3. Two different cannabis flower samples.



Figure 4. Six different cannabis flower samples.

Results and discussion

Calibration, accuracy, and precision

The main advantage of automated sample preparation is the reduction of errors and resources. The results show a robust process that can be easily implemented and maintained with minimal user intervention. Table 5 provides the retention time and linearity (R^2) for three independent calibration curves prepared using the PAL3 Series II RTC rail. All R^2 values were >0.99 , demonstrating excellent fit of prepared calibration standards. Table 6 displays the accuracy and precision data for three different calibration levels. All accuracy values fell between 91.5 to 118.6% and all precision values were under 8.7%.

Table 5. The retention time (RT) and linearity (R^2) for three independent calibration curves generated using the PAL3 Series II RTC rail. All calibration curves were fit with quadratic relationships, and 1/x weighting was used.

| Name | RT (min) | Cal 1 (R^2) | Cal 2 (R^2) | Cal 3 (R^2) |
|----------------------|----------|-----------------|-----------------|-----------------|
| alpha-Pinene | 13.963 | 0.9993 | 0.9980 | 0.9992 |
| Camphene | 14.710 | 0.9994 | 0.9984 | 0.9992 |
| Sabinene | 15.399 | 0.9996 | 0.9987 | 0.9989 |
| beta-Myrcene | 15.536 | 0.9989 | 0.9987 | 0.9991 |
| beta-Pinene | 15.665 | 0.9993 | 0.9984 | 0.9995 |
| alpha-Phellandrene | 16.449 | 0.9992 | 0.9989 | 0.9992 |
| delta-3-Carene | 16.496 | 0.9995 | 0.9984 | 0.9991 |
| alpha-Terpinene | 16.817 | 0.9982 | 0.9968 | 0.9978 |
| E-beta-Ocimene | 16.995 | 0.9996 | 0.9992 | 0.9987 |
| D-Limonene | 17.177 | 0.9995 | 0.9986 | 0.9988 |
| Z-beta-Ocimene | 17.490 | 0.9996 | 0.9991 | 0.9989 |
| Eucalyptol | 17.685 | 0.9998 | 0.9997 | 0.9992 |
| gamma-Terpinene | 18.072 | 0.9996 | 0.9989 | 0.9989 |
| Terpinolene | 19.053 | 0.9993 | 0.9984 | 0.9984 |
| Sabinene Hydrate | 19.484 | 0.9999 | 0.9998 | 0.9981 |
| Linalool | 19.779 | 1.0000 | 0.9998 | 0.9980 |
| [±]-Fenchone | 19.975 | 0.9998 | 0.9998 | 0.9989 |
| Endo-Fenchyl Alcohol | 20.446 | 0.9999 | 0.9997 | 0.9983 |
| Isopulegol | 20.802 | 0.9999 | 0.9998 | 0.9979 |
| [±/N]-Camphor | 21.088 | 0.9999 | 0.9998 | 0.9987 |
| Isoborneol | 21.186 | 0.9991 | 0.9999 | 0.9975 |
| Menthol | 21.207 | 0.9994 | 0.9996 | 0.9989 |
| [±]-Borneol | 21.330 | 0.9999 | 0.9997 | 0.9971 |
| alpha-Terpineol | 21.431 | 0.9999 | 0.9998 | 0.9981 |
| gamma-Terpineol | 21.431 | 0.9999 | 0.9998 | 0.9977 |
| Nerol | 21.729 | 0.9999 | 0.9999 | 0.9975 |
| Geraniol | 21.996 | 0.9999 | 0.9998 | 0.9972 |
| Pulegone | 22.147 | 0.9999 | 0.9998 | 0.9978 |
| Geranyl Acetate | 23.215 | 0.9999 | 0.9999 | 0.9985 |
| Farnesene | 23.935 | 0.9998 | 0.9989 | 0.9981 |
| alpha-Cedrene | 24.175 | 0.9999 | 0.9992 | 0.9987 |
| E-Caryophyllene | 24.227 | 0.9984 | 0.9990 | 0.9992 |
| alpha-Humulene | 24.721 | 0.9997 | 0.9996 | 0.9985 |
| Valencene | 25.106 | 0.9994 | 0.9995 | 0.9984 |
| Z-Nerolidol | 25.357 | 0.9999 | 0.9999 | 0.9991 |
| E-Nerolidol | 25.802 | 0.9999 | 0.9999 | 0.9991 |
| Guaiol | 27.097 | 0.9999 | 0.9999 | 0.9990 |
| Carophyllene Oxide | 27.460 | 0.9997 | 0.9998 | 0.9991 |
| Cedrol | 27.918 | 0.9999 | 1.0000 | 0.9992 |
| alpha-Bisabolol | 28.359 | 0.9999 | 0.9999 | 0.9990 |

Table 6. Accuracy as the average percent (n = 3) and precision as % RSD (n = 3) at calibration level 3.83, 30.64, and 245.10 µg/mL.

| Compound Name | Accuracy (average %, n = 3) | | | Precision (% RSD, n = 3) | | |
|----------------------|-----------------------------|-------------|--------------|--------------------------|-------------|--------------|
| | 3.83 µg/mL | 30.64 µg/mL | 245.10 µg/mL | 3.83 µg/mL | 30.64 µg/mL | 245.10 µg/mL |
| alpha-Pinene | 108.1 | 95.9 | 105.3 | 1.5 | 2.4 | 2.0 |
| Camphene | 108.0 | 95.9 | 104.8 | 1.4 | 2.2 | 1.7 |
| Sabinene | 108.5 | 95.4 | 104.5 | 1.0 | 1.6 | 1.7 |
| beta-Myrcene | 109.6 | 96.7 | 104.9 | 1.5 | 3.7 | 0.4 |
| beta-Pinene | 109.3 | 96.2 | 104.3 | 1.2 | 0.4 | 1.9 |
| alpha-Phellandrene | 105.8 | 97.8 | 102.0 | 1.4 | 4.1 | 4.0 |
| delta-3-Carene | 108.0 | 95.7 | 104.6 | 1.3 | 2.3 | 1.9 |
| alpha-Terpinene | 118.6 | 91.5 | 106.4 | 3.0 | 0.6 | 1.6 |
| E-beta-Ocimene | 111.2 | 95.2 | 104.0 | 3.8 | 1.4 | 1.2 |
| D-Limonene | 109.8 | 95.2 | 104.8 | 1.4 | 2.0 | 1.4 |
| Z-beta-Ocimene | 109.0 | 95.2 | 104.1 | 2.7 | 1.7 | 1.1 |
| Eucalyptol | 107.2 | 97.0 | 102.8 | 2.1 | 3.1 | 1.2 |
| gamma-Terpinene | 108.7 | 95.6 | 104.3 | 2.1 | 1.9 | 1.4 |
| Terpinolene | 112.5 | 94.3 | 105.0 | 2.4 | 2.0 | 1.3 |
| Sabinene Hydrate | 106.6 | 96.8 | 103.1 | 5.1 | 3.3 | 2.9 |
| Linalool | 105.9 | 96.5 | 103.2 | 5.1 | 3.6 | 3.0 |
| [±]-Fenchone | 106.9 | 96.7 | 103.0 | 3.6 | 2.5 | 1.5 |
| Endo-Fenchyl Alcohol | 107.7 | 96.8 | 103.1 | 1.5 | 2.1 | 2.8 |
| Isopulegol | 108.9 | 95.1 | 103.0 | 5.3 | 2.5 | 3.2 |
| [±/N]-Camphor | 107.3 | 97.0 | 102.9 | 5.3 | 3.6 | 1.9 |
| Isoborneol | 107.3 | 97.1 | 104.1 | 7.9 | 4.3 | 2.9 |
| Menthol | 107.5 | 97.7 | 103.4 | 4.0 | 3.7 | 1.2 |
| [±]-Borneol | 108.9 | 95.5 | 104.0 | 6.6 | 4.0 | 3.4 |
| alpha-Terpineol | 108.2 | 96.3 | 103.2 | 5.1 | 4.1 | 2.7 |
| gamma-Terpineol | 107.9 | 95.6 | 103.5 | 6.1 | 4.0 | 2.9 |
| Nerol | 105.9 | 98.0 | 103.2 | 8.7 | 5.4 | 3.5 |
| Geraniol | 111.2 | 95.1 | 103.4 | 7.6 | 5.4 | 3.4 |
| Pulegone | 108.6 | 96.1 | 103.4 | 6.3 | 4.2 | 2.8 |
| Geranyl Acetate | 106.7 | 97.1 | 102.9 | 3.2 | 3.5 | 2.3 |
| Farnesene | 109.1 | 98.6 | 104.1 | 7.5 | 4.4 | 2.4 |
| alpha-Cedrene | 108.6 | 94.6 | 103.5 | 4.8 | 3.1 | 1.8 |
| E-Caryophyllene | 107.0 | 97.0 | 104.9 | 3.7 | 1.4 | 0.9 |
| alpha-Humulene | 109.2 | 95.8 | 103.7 | 3.8 | 2.4 | 1.7 |
| Valencene | 108.2 | 95.5 | 104.3 | 4.3 | 2.2 | 1.4 |
| Z-Nerolidol | 107.7 | 96.2 | 102.1 | 3.0 | 2.7 | 1.8 |
| E-Nerolidol | 106.3 | 96.6 | 102.1 | 3.5 | 2.5 | 1.8 |
| Guaiol | 108.0 | 96.3 | 102.1 | 3.7 | 2.0 | 2.0 |
| Carophyllene Oxide | 108.8 | 95.4 | 102.7 | 1.2 | 1.6 | 1.4 |
| Cedrol | 106.1 | 96.4 | 101.7 | 3.6 | 2.0 | 2.0 |
| alpha-Bisabolol | 106.7 | 96.3 | 102.5 | 4.4 | 2.3 | 1.7 |

Analysis of various cannabis samples

All limits of detection (LOD) and limits of quantitation (LOQ) fell below the lowest calibration point (3.83 µg/mL). Thus, the reporting limit for the current study was defined as any value greater than 3.83 µg/mL. The terpene concentrations found in the different cannabis samples are presented in Table 7.

Each sample shows a unique terpene profile. Concentration of detected terpenes in four replicates are presented for sample 5 (Figure 5) and sample 6 (Figure 6). The % RSD for detected terpenes ranged between 0.69 to 7.92% in sample 5 and 3.47 to 10.16% in sample 6.

Table 7. Concentration (µg/mL) of terpenes found in 6 different cannabis flower samples. Samples 5 and 6 were run in quadruplicate (n = 4) to determine the standard deviation (SD).

| Compound Name | Sample 1 (µg/mL) | Sample 2 (µg/mL) | Sample 3 (µg/mL) | Sample 4 (µg/mL) | Sample 5 (µg/mL ±SD, n = 4) | Sample 6 (µg/mL ±SD, n = 4) |
|----------------------|------------------|------------------|------------------|------------------|-----------------------------|-----------------------------|
| alpha-Pinene | N/A | 4.20 | N/A | N/A | N/A | N/A |
| D-Limonene | N/A | N/A | N/A | N/A | 4.08 ±0.12 | N/A |
| Eucalyptol | 8.02 | N/A | N/A | N/A | N/A | N/A |
| Linalool | 19.26 | 4.45 | 9.78 | 4.23 | 10.16 ±0.44 | N/A |
| Endo-Fenchyl Alcohol | 7.46 | N/A | 11.26 | N/A | 8.22 ±0.06 | N/A |
| [±]-Borneol | N/A | N/A | 7.72 | N/A | N/A | N/A |
| alpha-Terpineol | 18.39 | N/A | 35.21 | 7.57 | 9.26 ±0.45 | 4.02 ±0.25 |
| gamma-Terpineol | 15.34 | N/A | 24.71 | 6.90 | 8.18 ±0.47 | N/A |
| Farnesene | 139.52 | 20.34 | 98.73 | 130.25 | 76.83 ±6.08 | 31.24 ±3.17 |
| E-Caryophyllene | 95.40 | 21.35 | 57.19 | 63.06 | 159.99 ±4.61 | 21.12 ±0.73 |
| alpha-Humulene | 38.52 | 7.87 | 20.82 | 17.61 | 40.69 ±1.41 | 6.33 ±0.50 |
| E-Nerolidol | 13.92 | N/A | 14.50 | 5.38 | 7.18 ±0.38 | N/A |
| Guaiol | N/A | N/A | 9.04 | N/A | N/A | N/A |
| Carophyllene Oxide | 26.59 | 20.78 | 50.75 | 44.05 | 20.97 ±1.08 | 8.47 ±0.40 |
| alpha-Bisabolol | 16.82 | 29.43 | 48.34 | 21.02 | 117.33 ±5.29 | 23.23 ±0.94 |

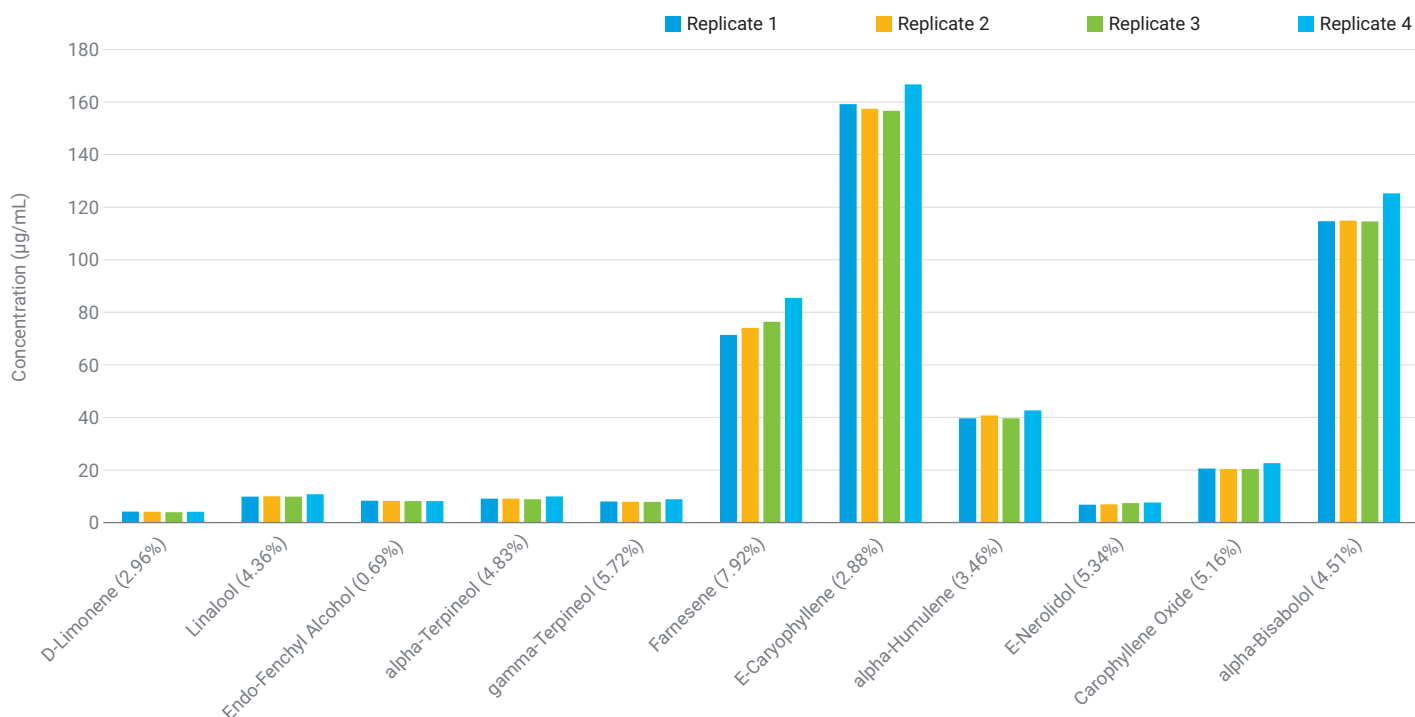


Figure 5. Terpenes concentration (µg/mL) in sample 5 (n = 4). The % RSD among replicates is shown in parenthesis.

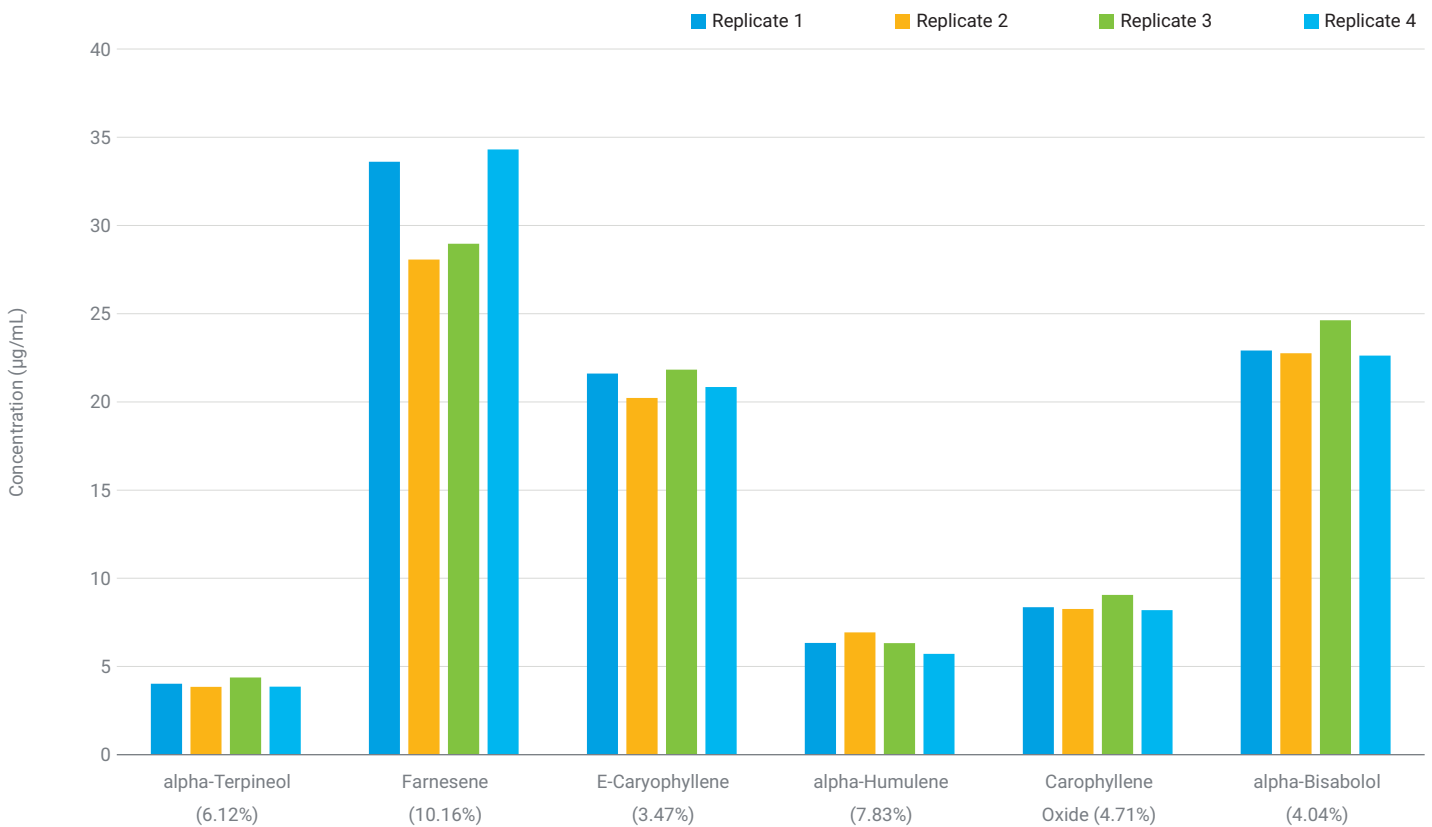


Figure 6. Terpenes concentration (µg/mL) in sample 6 (n = 4). The % RSD among replicates is shown in parenthesis.

Conclusion

The presented application demonstrates that the Agilent PAL3 Series II RTC sampler can accurately and precisely conduct the automated sample preparation for the analysis of terpenes in cannabis flower. The use of automated sample preparation helps reduce user errors and the amount of sample and solvent needed for analysis. Coupled with an Agilent Intuvo 9000/5977C GC/MS system, PAL3 Series II RTC automation can save time and reduce energy consumption, leading to a greener high-throughput laboratory.

Reference

- Hollis, J. S.; Harper, T.; Macherone, A. Terpenes Analysis in Cannabis Products by Liquid Injection using the Agilent Intuvo 9000/5977B GC/MS System, *Agilent Technologies application note*, publication number 5994-2032EN.

Agilent products and solutions are intended to be used for cannabis quality control and safety testing in laboratories where such use is permitted under state and country law.

www.agilent.com

DE40874749

This information is subject to change without notice.

© Agilent Technologies, Inc. 2023
 Printed in the USA, June 7, 2023
 5994-6007EN