

# Terpenes and residual solvents in hemp products by GC-FID



Complete CANNABIS TESTING SOLUTIONS for the Canadian market from one trusted source

Authors: CCS and GC Center of Excellence (CoE) Application Scientists, Thermo Fisher Scientific

Keywords: Hemp, terpenes, residual solvents

## Introduction

The production of hemp extracts and concentrates requires the use of solvents such as ethanol, butane or propane. The presence of these solvents in final consumer products must be monitored for safety and quality purposes.

Terpenes are the compounds that provide the unique aroma of hemp and hemp products and are thought to impact the user's experience. Although not typically regulated, terpene profiles can provide valuable information for product quality.

This protocol provides a starting point for the development of a method for detection and quantitation of terpenes and residual solvents in hemp and hemp products for GC analysis with headspace sampling.

## Important notes

- This application brief provides starting conditions for residual solvents and terpenes on the Thermo Scientific™ TraceGOLD™ TG-624SiIMS GC column, allowing for users to optimize their chromatography for desired target analytes and limits, depending on the local rules and regulations



- For plant material and many oils, concentrates and extracts, a small aliquot (for example, 10 mg or 10  $\mu$ L) may be placed directly into the headspace vial. If needed, samples may be dissolved in dimethylacetamide (DMA) or another appropriate solvent before analysis.
- The high thermal stability of the TraceGOLD TG-624SiIMS GC columns allow for consistent chromatography of late-eluting terpenes

## Materials required

- GC-FID instrument such as Thermo Scientific™ TRACE™ 1310 Gas Chromatograph with headspace autosampler
- TraceGOLD TG-624SiIMS GC Column, 60 m  $\times$  0.32 mm ID  $\times$  1.8  $\mu$ m (P/N 26059-3410)

## Protocol

1. Analyze samples using the following GC conditions:

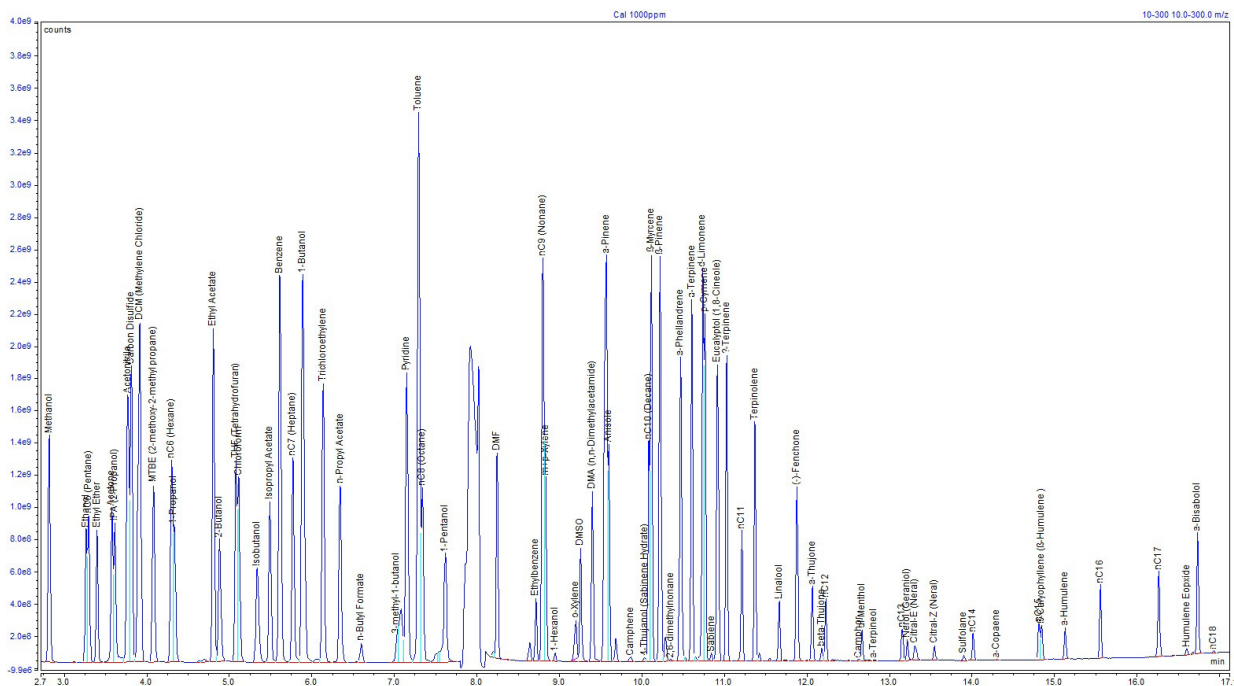
Headspace parameters	Incubation: 80 °C for 2.5 min
	Syringe temp: 90 °C
	Injection volume: 0.7 mL, Filling vol.: 0.8 mL
	Purge: N <sub>2</sub> , 5 s pre-injection, 60 s post-injection
Inlet module mode	SSL, Split with cyclosplitter inlet liner
Split ratio	10:1
Inlet temperature	225 °C
Carrier gas, mode, flow	Helium, 3.25 mL/min constant flow
Oven program	40 °C initial, hold 0.2 min
	17.5 °C/min to 90 °C, hold 1.5 min
	17.5 °C/min to 310 °C, hold 1.37 min
	Total run time of 18.5 min
FID temperature	310 °C

- Most residual solvents elute in the first half of the chromatogram. If these are the only compounds of interest the run can be shortened.
- Terpenes begin to elute at approximately 10 minutes in this method. A higher initial temperature can be used for analysis of terpenes only.

## Residual solvents and terpenes RTs

Compound	Approximate retention time (min)
Methanol	2.82
Ethanol	3.27
nC5 (Pentane)	3.30
Acetone	3.58
IPA (2-Propanol)	3.62
Acetonitrile	3.78
Carbon Disulfide	3.82
DCM (Methylene Chloride)	3.92
MTBE (2-methoxy-2-methyl propane)	4.09
nC6 (Hexane)	4.31
MEK (2-Butanone)	4.81
2-Butanol	4.89
THF (Tetrahydrofuran)	5.09
Chloroform	5.12
Benzene	5.62
nC7 (Heptane)	5.78
1-Butanol	5.89
Trichloroethylene	6.14
n-Butyl Formate	6.60
Pyridine	7.16
Toluene	7.30
Ethylbenzene	8.72
m+p-Xylene	8.84
o-Xylene	9.21
DMSO	9.26
DMA (n,n-Dimethylacetamide)	9.40
a-Pinene	9.57
Camphene	9.87
β-Pinene	10.23

Compound	Approximate retention time (min)
a-Phellandrene	10.47
a-Terpinene	10.61
d-Limonene	10.74
p-Cymene	10.77
Sabien	10.85
Eucalyptol (1,8-Cineole)	10.92
Terpinene	11.03
nC11	11.22
Terpinolene	11.37
Linalool	11.67
(-)-Fenchone	11.88
a-Thujone	12.07
beta-Thujone	12.18
nC12	12.23
dl-Menthol	12.67
nC13	13.16
Nerol (Geraniol)	13.22
Citral-Z (Neral)	13.55
Sulfolane	13.90
nC14	14.01
nC15	14.81
β-Caryophyllene (β-Humulene )	14.84
a-Humulene	15.13
nC16	15.56
nC17	16.27
Humulene Epoxide	16.61
a-Bisabolol	16.74
nC18	16.93



Related products

Description	Part number
TraceGOLD TG-624SiIMS column, 60 m × 0.32 mm ID × 1.8 μm	26059-3410
Amber headspace glass vial with crimp cap, 10 mL	10-CV(A)
Headspace glass vial with blue magnetic crimp cap with seal	20-MCBC-ST3



Offering one of the broadest portfolios of cannabis testing equipment for the Canadian cannabis market, Thermo Fisher Scientific is your one-stop resource for customized solutions, proven workflows and support. Whether setting up your first laboratory or scaling existing operations, we offer a complete range of instrument workflows to help you comply with Canadian cannabis testing regulations.

Find out more at [thermofisher.com/cannabistesting](https://thermofisher.com/cannabistesting)

Current versions of product instructions are available at [thermofisher.com/chromexpert](https://thermofisher.com/chromexpert)

Find more information about GC columns at [thermofisher.com/gccolumns](https://thermofisher.com/gccolumns)

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. Thermo Fisher Scientific does not support, encourage or promote the use of its products or services in connection with any illegal use, cultivation or trade of cannabis or cannabis products. Thermo Fisher Scientific products are intended to be used only in compliance with all applicable laws in a manner that promotes public safety and/or in connection with any lawful and approved scientific or medical research activities. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all locations. Please consult your local sales representatives for details. **AB22001-EN 0520M**