

GC-IRMS

GC-IRMS: Benefits of Static Headspace Sampling for carbon isotope analysis of methanol and ethanol in water matrix; applicability to wine and spirits ethanol $\delta^{13}\text{C}$ determination

Authors

María de Castro¹, Ana I. Cabañero²,
Vanessa Peiro³, Mario Tuthorn¹

¹Thermo Fisher Scientific

²Laboratorio Arbitral Agroalimentario,
Ministerio de Agricultura, Pesca y
Alimentación (MAPA), Spain

³SIdI, Universidad Autónoma de Madrid
(UAM), Spain

Keywords

VOC, Alcohol, Wine, Food, SHS
sampling, IRMS, isotope fingerprints

Introduction

Analyzing alcohols in water matrix using gas chromatography (GC) is challenging due to the large aqueous component content in the sample matrix. Water in the sample shortens the lifespan of GC columns and requires frequent maintenance of the injector and column. This can complicate the analysis of isotope signatures in wine and spirits by GC-C-IRMS, for example, which is crucial for verifying product authenticity and preventing fraud. In this case, the methodology focuses on $\delta^{13}\text{C}$ measurements of ethanol (EtOH), a useful indicator for detecting wine and spirits adulterations and determining its origin.^{1,2}

Measuring $\delta^{13}\text{C}$ content can detect and quantify the addition of C4 plant sugars (such as sugar cane or corn isoglucose) to grape-derived products. The current official analytical procedures face challenges such as multiple steps required to extract EtOH from the wine matrix (e.g., distillation) and technical difficulties in ensuring the collected EtOH is free from isotope fractionation effects.

An optimized method for carbon isotope analysis of volatile alcohols in water matrix by GC-IRMS aims to streamline the process by eliminating the need for prior alcohol isolation, thus simplifying sample preparation and speeding up analysis. This technical note aims to demonstrate that Static Headspace Sampling (SHS) injection of methanol (MeOH) and EtOH in water matrix via a split/splitless injector simplifies not only sample preparation, but reduces column contamination, minimizes system maintenance, and produces robust and high-quality isotopic data. We report optimized methodology based on EtOH and MeOH standards, and EtOH $\delta^{13}\text{C}$ isotopic data for wine and spirits samples.



Analytical setup

All measurements are performed using a Thermo Scientific™ GC IsoLink™ II IRMS System with the Thermo Scientific™ TriPlus™ RSH™ Series Autosampler, equipped with a Static Headspace Sampling (SHS) injection option (Figure 1).

For the SHS technique, the samples are placed in tightly closed vials. Each vial is heated, and the volatile compounds are transferred from the liquid sample into the gaseous phase above it (headspace) until a condition of thermodynamic equilibrium is reached. Afterwards, an aliquot of headspace is withdrawn and injected into the gas chromatograph. The trend of a volatile compound to transfer into the headspace is expressed by the partition coefficient *K*. The coefficient depends on the compound solubility in the matrix, and it is strongly affected by the temperature and the composition of the matrix itself. The analytical times are optimized according to the set parameters, such as equilibration time and analytical time, to fully utilize the instrument capability.

EtOH and MeOH isotopically certified standards from the Indiana University were used for:

1. Developing chromatography method
2. Testing repeatability of the results



Figure 1. TriPlus RSH Series Autosampler equipped with a Static Headspace Sampling (SHS) injection option.

3. Confirming that the developed methodology is delivering accurate isotopic data
4. Defining LOD for each compound and evaluating possible linearity effects

The standards stock solutions were prepared as outlined in Table 1.

Table 1. Dilution of alcohol standards for carbon isotope analysis using GC-C-IRMS with Static Headspace Sampling injection

| Name (Manufacturer) | Components | Original solution | Stock solution ¹ | Standards preparation in salty water ² |
|-------------------------------------|-----------------------------------|---|---|---|
| MeOH (Indiana University) | MeOH | ≥ 99.9 wt. % | 20 mL H ₂ O Milli-Q + 76 µL MeOH = 3000 ppm stock solution | 20 mL vial: 9.5 mL Milli-Q H ₂ O with NaCl 5.1 M + 500 µL std solution = approx. 150 ppm MeOH |
| EtOH #3 *(Indiana University) | EtOH distillate from vodka | 82 wt. % (rest water) | 20 mL H ₂ O Milli-Q + 76 µL EtOH = 3000 ppm stock solution | 20 mL vial: 9.5 mL Milli-Q H ₂ O with NaCl 5.1 M + 500 µL std solution = approx. 150 ppm EtOH #3 |
| EtOH #4 *(Indiana University) | EtOH distillate from rum | 80.7 wt. % (rest water) | 20 mL H ₂ O Milli-Q + 76 µL EtOH = 3000 ppm stock solution | 20 mL vial: 9.5 mL Milli-Q H ₂ O with NaCl 5.1 M + 500 µL std solution = approx. 150 ppm EtOH #4 |
| MeOH + EtOH #3 (Indiana University) | MeOH + EtOH distillate from vodka | ≥ 99.9 wt. % // 82 wt. % (rest water) | na | 20 mL vial: 9 mL Milli-Q H ₂ O with NaCl 5.1 M + 500 µL stock MeOH + 500 µL stock EtOH #3 = approx. 150 ppm each |
| MeOH + EtOH #4 (Indiana University) | MeOH + EtOH distillate from rum | ≥ 99.9 wt. % // 80.7 wt. % (rest water) | na | 20 mL vial: 9 mL Milli-Q H ₂ O with NaCl 5.1 M + 500 µL stock MeOH + 500 µL stock EtOH #4 = approx. 150 ppm each |

¹MeOH density is 0.792Kg/L and EtOH density 0.790Kg/L, which is considered for concentration calculations of the standards solutions. Ethanol #3 is an azeotropic distillate from vodka (C3) with 82 wt. % (87.32 vol. %), and Ethanol #4 is a rum (C4) azeotropic distillate with 80.7 wt. %, but this is not considered for concentration calculations of the solutions.

²To lower partitioning coefficient of target compounds, it is recommended to add approximately 3 g of NaCl to every vial before closure. This promotes higher concentration of VOCs in the gas phase and helps avoiding fractionation due to lighter isotopic composition in the gas phase. To facilitate the standards and samples preparations, a solution of Milli Q water with NaCl 5.1M was prepared by adding 300 grams of NaCl to one liter of Milli Q water, which avoids the manual addition of NaCl to each vial.

*NOTE: There are two new and purer EtOH standards available from Indiana University: EtOH#1 from C3 plant (99.96 vol %) and Ethanol #2 from C4 plant (99.11 vol. %).

For the evaluation of the possible linearity effects, concentration gradient solutions were prepared using the 3000 ppm stock solution.

Detailed analytical setup for the TriPlus RSH Series Autosampler with the Static Headspace Sampling option, the Thermo Scientific™ iConnect™ Split/Splitless (SSL) Injector

Module and the TRACE™ Series GC is listed in Table 2. The GC IsoLink II IRMS System operation is driven by Thermo Scientific™ Qtegra™ ISDS Software.

Conditioning of the combustion reactor in the GC IsoLink II conversion interface was performed as outlined in Table 3.

Table 2. Autosampler and gas chromatographer analytical setup for the isotopic analysis of MeOH and EtOH

| Triplus RSH Series autosampler parameters | | iConnect SSL Injector Module parameters | |
|---|------------------|--|--------------------|
| Thermo Scientific™ Fixed Needle Autosampler Syringe, Gas-tight, Headspace, 2.5 mL, 23 G, 65 mm (PN 365Q2131/PN 365L2321-SM) | | Thermo Scientific™ LinerGOLD™ GC Liners, Direct Straight Liner (PN 453A1335-UI) | |
| Injection type | Static Headspace | Injection temperature | 150 °C |
| Sample draw | 0.5 mL | Inlet module and mode | SSL, split |
| Sampling depth mode | Standard | Split flow | 24 mL/min |
| Agitator temperature | 65 °C | Split ratio | 20:1 |
| Incubation time | 15 min | Septum purge flow | 3 mL/min, constant |
| Agitation speed | 250 rpm | Carrier gas, flow | Helium, 1.2 mL/min |
| Agitator on | 10 sec | TRACE Series GC parameters | |
| Agitator off | 2 sec | Thermo Scientific™ TraceGOLD™ TR-WAX GC Column, 60 m, 0.25 mm, 0.25 µm (PN 260W154P) | |
| Syringe temperature | 70 °C | Temperature | 40 °C |
| Fill strokes volume | n.a. | Hold time | 3.5 min |
| Fill strokes count | 0 | Rate | 15 °C/m |
| Filling delay | 3 sec | Temperature 2 | 95 °C |
| Pre-injection syringe flush | Enabled | Hold time | 2 min |
| Post-injection syringe flush | 60 s | Rate 2 | 25 °C/min |
| Filling speed | 10 mL/min | Temperature 3 | 220 °C |
| Injection speed | 30 mL/min | Hold time | 4 min |
| Injection depth | 45 mm | GC run time | 18.2 min |
| Penetration speed | 25 mm/s | Prep run timeout | 120 min |
| Pre-injection delay | 1 s | | |
| Post-injection delay | 3 s | | |
| Needle speed in vial | 20 mm/s | | |
| Synchro type | Normal | | |

Table 3. Conditioning parameters for the combustion reactor of the GC IsoLink II Conversion Unit

| Reactor conditioning type | Setup |
|--|--|
| Extended – used at the start of a sequence | 45 min O ₂ + 90 min Backflush + 0.1 min purge |
| Seed – used with each sample run | 0.2-0.3 min O ₂ + 0.2 min Backflush + 0.1 min purge |

For optimization of the SHS injection of MeOH and EtOH via split/splitless injector the following analytical steps were followed:

1. Individual alcohols standards injections were performed for developing the chromatography method, testing repeatability and accuracy of the results
2. A certified standards mix (MeOH + EtOH #3, from Indiana University) was prepared to evaluate its chromatographic separation, as well as to ensure that the precision and accuracy of the data were not compromised by injecting both alcohols together, since you could find both of them at the same time in real samples

3. Concentration gradients were analyzed for the evaluation of possible linearity effects using isotopically certified standards (Indiana University)

Results

For the analysis, 9.5 mL of Milli-Q H₂O with NaCl 5.1 M and 500 µL of stock solution were transferred into 20 mL, 18 mm screw top, headspace vials (PN 6ASV20-1) with 18mm magnetic screw caps (PN 6PMSC18-STH, 8 mm hole). Chromatographic separation of the target compounds is shown in Figure 2.

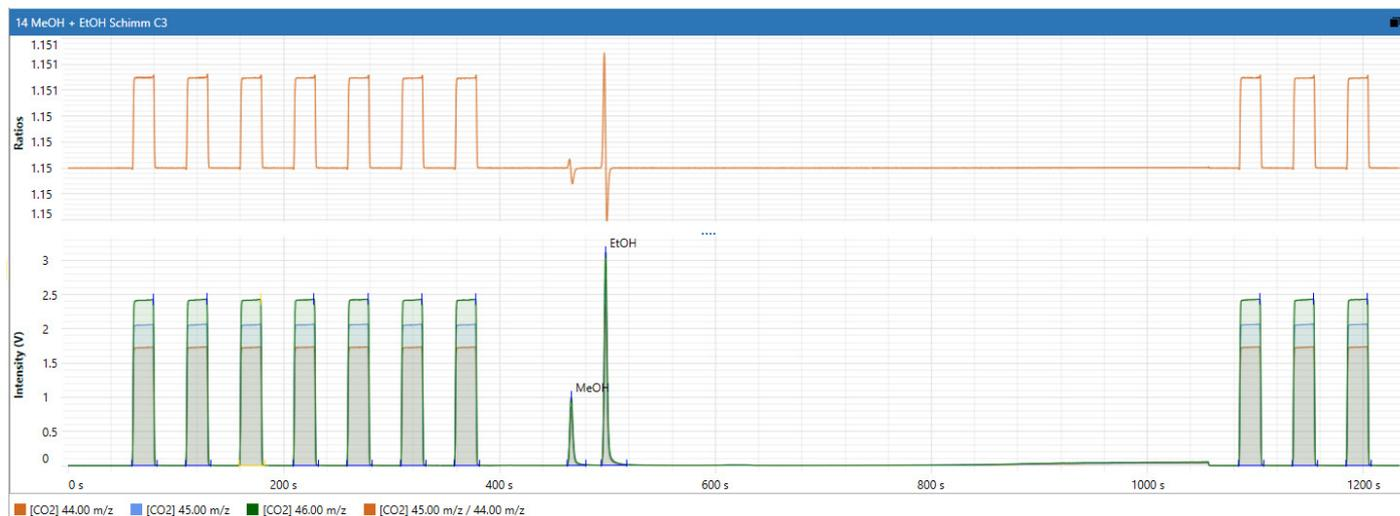


Figure 2. GC-IRMS chromatogram of the MeOH and EtOH #3 certified standards mix; 150 ng/uL (150ppm) each in H₂O

Data evaluation was performed using Qtegra ISDS Software tools for data normalization, where first isotopically certified standards injections were treated as Compound Specific Isotope Analysis (CSIA) δ Standard and the following 10

repetitions were treated as unknown samples (see Sample list in Figure 3). The resulting in-software evaluation shows high correlation between the measured and known isotope ratios values (Figure 4).

| Sample List | Label | Status | Comment | Evaluate | Sample Type | Reference | Vial | Inject Volume [µl] | Inject |
|-------------|-----------------------|--------|-----------------------------|----------|-----------------------|--------------------|------|--------------------|--------|
| 1 | Blank | ● | 45mins O2 + 2hs BF_18s seed | ✓ | Unknown | | 164 | 500 | ✓ |
| 2 | MeOH + EtOH Schimm C3 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Delta Standard (CSIA) | Alcohols Schimm #3 | 171 | 500 | ✓ |
| 3 | EtOH Schimm C4 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Delta Standard (CSIA) | EtOH Schimm #4 | 172 | 500 | ✓ |
| 4 | MeOH + EtOH Schimm C3 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 173 | 500 | ✓ |
| 5 | EtOH Schimm C4 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 174 | 500 | ✓ |
| 6 | MeOH + EtOH Schimm C3 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 175 | 500 | ✓ |
| 7 | EtOH Schimm C4 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 176 | 500 | ✓ |
| 8 | MeOH + EtOH Schimm C3 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 177 | 500 | ✓ |
| 9 | EtOH Schimm C4 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 178 | 500 | ✓ |
| 10 | MeOH + EtOH Schimm C3 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 179 | 500 | ✓ |
| 11 | EtOH Schimm C4 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 180 | 500 | ✓ |
| 12 | MeOH + EtOH Schimm C3 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 181 | 500 | ✓ |
| 13 | EtOH Schimm C4 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 182 | 500 | ✓ |
| 14 | MeOH + EtOH Schimm C3 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 183 | 500 | ✓ |
| 15 | EtOH Schimm C4 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 184 | 500 | ✓ |
| 16 | MeOH + EtOH Schimm C3 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 185 | 500 | ✓ |
| 17 | EtOH Schimm C4 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 186 | 500 | ✓ |
| 18 | MeOH + EtOH Schimm C3 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 187 | 500 | ✓ |
| 19 | EtOH Schimm C4 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 188 | 500 | ✓ |
| 20 | MeOH + EtOH Schimm C3 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 189 | 500 | ✓ |
| 21 | EtOH Schimm C4 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 190 | 500 | ✓ |
| 22 | MeOH + EtOH Schimm C3 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 191 | 500 | ✓ |
| 23 | EtOH Schimm C4 | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 192 | 500 | ✓ |
| 24 | Blank | ● | 5ml in 10ml vial_1.5g NaCl | ✓ | Unknown | | 165 | 500 | ✓ |

Figure 3. Qtegra ISDS Software sample list for in-software data normalization using a CSIA standard

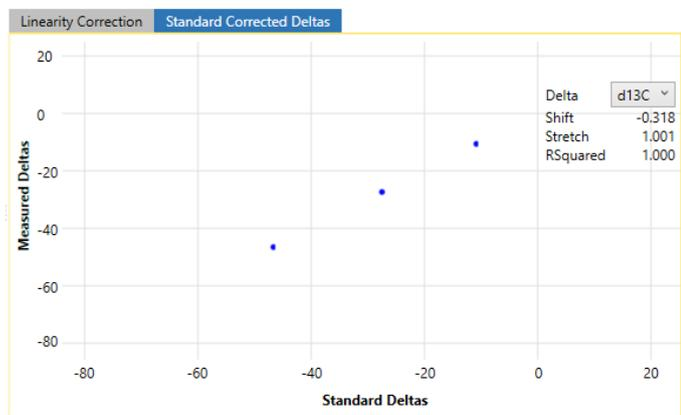


Figure 4. Qtegra ISDS Software data evaluation of measured isotope values vs. known isotope values of a CSIA standard

The certified standards isotope data reported in Table 4 demonstrate excellent measurement precision and accuracy for all compounds, as well as good RSD for total areas.

The Figure 5 shows data from three different analysis sequences done during the method development process, showing excellent intra and inter-sequence repeatability, and confirming that the developed methodology is delivering accurate isotopic data.

Table 4. Isotope data for the MeOH and EtOH standards; 150 ng/μL (150 ppm) in H₂O acquired using an SHS injection option

| Component | δ ¹³ C (‰) | | | | Total Area (V/s) | | |
|-----------|---------------------------------|--------------------|-----------|-------------------------|------------------|-----------|---------|
| | Certified δ ¹³ C (‰) | Average (‰) (n=10) | SD (n=10) | Offset vs Certified (‰) | Average (n=10) | SD (n=10) | RSD (%) |
| MeOH | -46.77 | -46.75 | 0.10 | 0.02 | 6.76 | 0.20 | 2.97 |
| EtOH #3 | -27.53 | -27.43 | 0.07 | 0.10 | 15.63 | 0.39 | 2.48 |
| EtOH #4 | -10.98 | -11.05 | 0.08 | -0.07 | 13.68 | 0.41 | 3.01 |

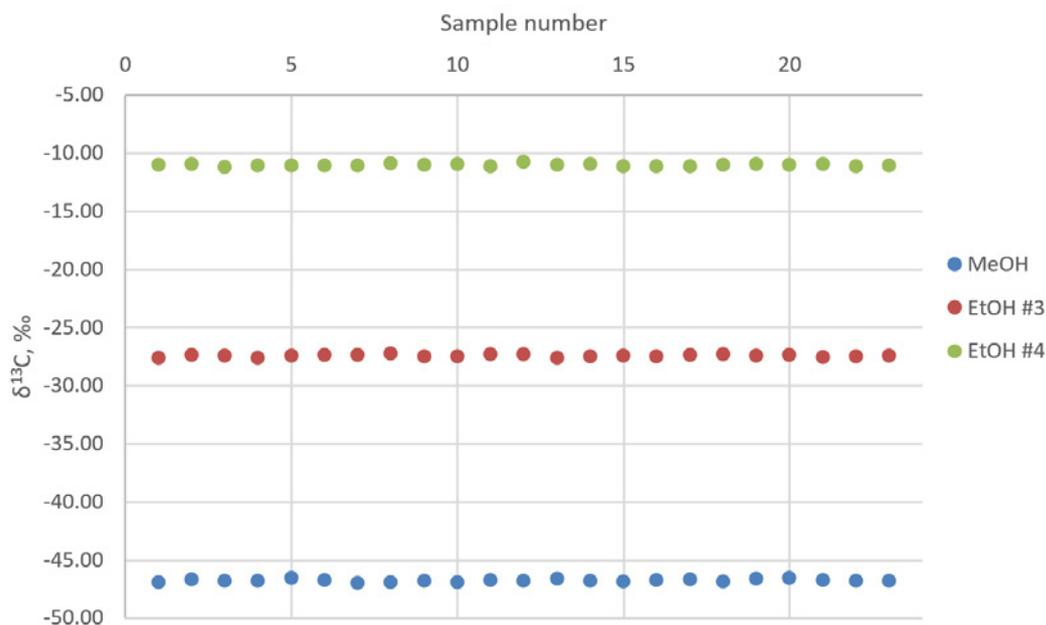


Figure 5. Isotope data acquired in different sequences for the MeOH and EtOH certified standards; 150 ng/μL (150 ppm) in H₂O using an SHS injection option

Finally, concentration gradients of the MeOH + EtOH #4 certified standard mix were analyzed with results showing that the isotope ratio of MeOH and EtOH are independent of the

amount of material analyzed within the range from 120 ppm to 1500 ppm (Table 5).

Table 5. Isotope ratio data of MeOH and EtOH #4 mix in the concentration range from 120 ppm to 1500 ppm

| Compound | MeOH | | | | | EtOH #4 | | | | |
|-------------------------------------|---------------|----------|---------------------------|----------|-----------------|---------------|----------|---------------------------|----------|-----------------|
| Certified $\delta^{13}\text{C}$ (‰) | -46.77 | | | | | -10.98 | | | | |
| ng/ μL (ppm) | Ampl. 44 (mV) | | $\delta^{13}\text{C}$ (‰) | | | Ampl. 44 (mV) | | $\delta^{13}\text{C}$ (‰) | | |
| | Avg. (n=3) | SD (n=3) | Avg. (n=3) | SD (n=3) | Offset vs Cert. | Avg. (n=3) | SD (n=3) | Avg. (n=3) | SD (n=3) | Offset vs Cert. |
| 120 | 0.32 | 0.02 | -46.95 | 0.03 | -0.18 | 1.03 | 0.06 | -10.78 | 0.28 | 0.20 |
| 240 | 0.66 | 0.01 | -46.75 | 0.05 | 0.02 | 2.08 | 0.02 | -11.19 | 0.13 | -0.21 |
| 480 | 1.34 | 0.08 | -46.51 | 0.14 | 0.26 | 4.15 | 0.23 | -10.68 | 0.21 | 0.30 |
| 1000 | 2.58 | 0.18 | -46.66 | 0.13 | 0.11 | 7.71 | 0.47 | -10.75 | 0.07 | 0.23 |
| 1500 | 3.88 | 0.13 | -46.57 | 0.12 | 0.20 | 10.82 | 0.26 | -10.79 | 0.01 | 0.19 |

Analysis of EtOH in wine and spirits samples

The Laboratorio Arbitral Agroalimentario – MAPA, Spain, supplied an in-house wine standard (Quality Control, QC) and wine and spirits samples that had been previously analyzed using the OIV official method (ethanol collection from wine with Cadiot columns for EA-IRMS analysis). These were provided along with each sample distillate for comparison with the SHS GC-IRMS methodology.

Some of the wine samples may still contain active bacteria. To prevent potential bacterial activity, two valid procedures can be used: filtration (0.22 μm) or sterilization. For this study, sterilization with benzoic acid was chosen. To determine the required amount and effectiveness of benzoic acid, two vials of each wine type (white and red) were prepared. One vial, without benzoic acid, was analyzed immediately after preparation. The other vial, containing approximately 25 mg of benzoic acid (a spatula tip), was stored at room temperature for 24 hours before analysis. The results (peak size and $\delta^{13}\text{C}$) were identical, demonstrating 100% efficacy.

Note: the total sample volume can be filtered when transferring it to the 2 mL storage vial, or the entire sample volume can be sterilized directly in the storage vial. This eliminates the need to add benzoic acid to each individual vial.

To simplify the preparation of standards and samples, a solution of Milli-Q water with 5.1 M NaCl was prepared by adding 300 grams of NaCl to one liter of Milli-Q water. This avoids the need to manually add NaCl to each vial. Stock solutions of the isotopically certified standards and samples distillates were prepared as shown in Table 6, with concentrations similar to those expected in the samples.

According to the principle of identical treatment, benzoic acid should be added to the standards and samples distillates vials as well, just as it is added to the wine and spirits samples. However, a test was conducted to evaluate any possible effects of benzoic acid addition by analyzing standards prepared with and without benzoic acid. The results showed no effect, therefore it is not necessary to add benzoic acid to the standards and samples distillates vials.

Table 6. Stock solutions of the isotopically certified standards and samples distillates to be analysed along with the wine and spirits samples

| Name (Manufacturer) | Components | Original solution | Stock solution in H_2O with NaCl 5.1 M |
|-------------------------------------|-----------------------------------|--|--|
| EtOH #4 (Indiana University) | EtOH distillate from rum | 80.7 wt. % (rest water) | 2 mL vial: 360 μL EtOH + 1640 μL Milli-Q H_2O = 150000 ppm (15 vol. %) |
| MeOH + EtOH #3 (Indiana University) | MeOH + EtOH distillate from vodka | ≥ 99.9 % // 82 wt. % (rest water) | 2 mL vial: MeOH 300 μL + EtOH 360 μL + 1340 Milli-Q H_2O = 150000 ppm (15 vol. %) each |
| Samples distillates | EtOH distillate from samples | 92 wt. % (rest water) | 2 mL vial: 360 μL EtOH + 1640 μL Milli-Q H_2O = 150000 ppm (15 vol. %) |

For the analysis, 10 mL of Milli-Q H₂O with 5.1 M NaCl were transferred into 20 mL, 18 mm screw top, headspace vials with 18 mm magnetic screw caps. Approximately 25 mg of benzoic acid was added to each vial that will be containing wine or spirits samples. Then, 50 µL of wine or distillate, and stock standard solutions were added to the vials. For the spirits

samples, a volume between 25-40 µL was added to the vial (depending on the EtOH vol. %).

Figure 6 shows clean separation of EtOH from other components in a wine sample analyzed by SHS GC-IRMS method.

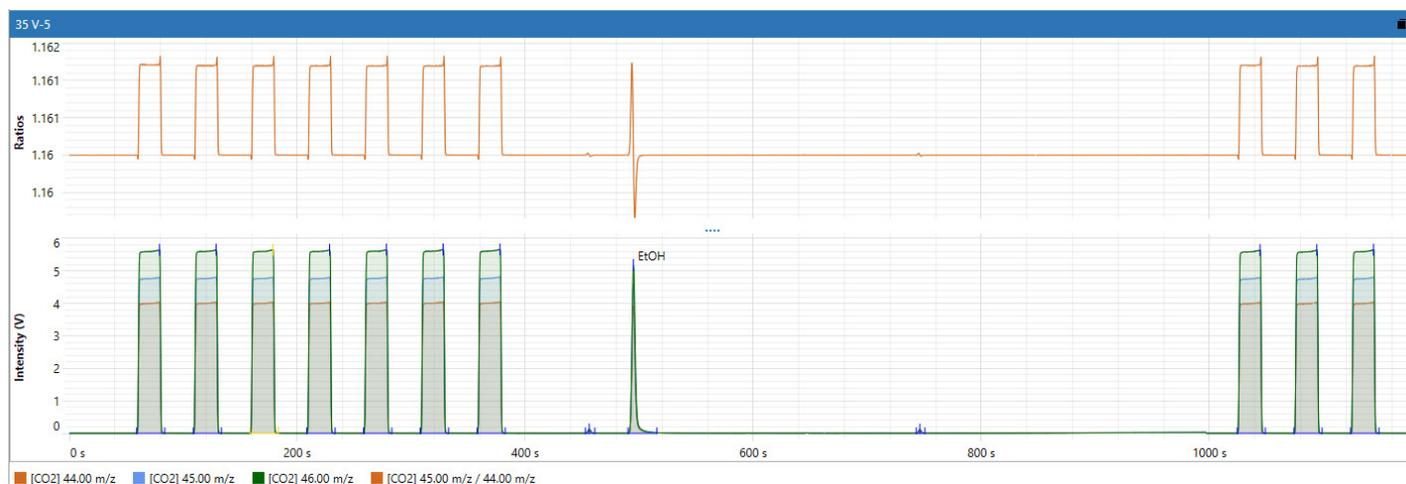


Figure 6. GC-IRMS chromatogram of the EtOH in a wine sample using SHS methodology

Table 7 shows the results obtained by SHS GC-IRMS and comparison with the data obtained from the Laboratorio Arbitral Agroalimentario – MAPA, Spain, for the samples

distillates analysis by EA-IRMS using the official method. We report excellent correlation of the SHS GC-IRMS and EA-IRMS data.

Table 7. Carbon isotope data of wine and spirits samples and samples distillates acquired by SHS GC-IRMS methodology and comparison to EA-IRMS data

| Sample | Sample | Volume added to vial (µL) | Average SHS GC-IRMS $\delta^{13}\text{C}$ (‰) (n=3) | SD (‰) (n=3) | EA-IRMS MAPA $\delta^{13}\text{C}$ (‰) | Offset GC-IRMS vs EA-IRMS (‰) |
|-----------------------|-----------------------------------|---------------------------|---|--------------|--|-------------------------------|
| Orujo (~40 vol. %) | Distillate stock (EtOH 15 vol. %) | 50 | -26.52 | 0.08 | -26.56 | -0.04 |
| | Raw sample | 25 | -26.51 | 0.10 | | -0.05 |
| Vermut (~15 vol. %) | Distillate stock (EtOH 15 vol. %) | 50 | -26.69 | 0.08 | -26.70 | -0.01 |
| | Raw sample | 40 | -26.82 | 0.11 | | 0.12 |
| Brandy (~40 vol. %) | Distillate stock (EtOH 15 vol. %) | 50 | -26.48 | 0.08 | -26.26 | 0.22 |
| | Raw sample | 25 | -26.54 | 0.10 | | 0.27 |
| Red wine (~14 vol. %) | Distillate stock (EtOH 15 vol. %) | 50 | -25.69 | 0.08 | -25.62 | 0.07 |
| | Raw sample | 50 | -25.71 | 0.07 | | 0.09 |
| Wine QC (~14 vol. %) | Distillate stock (EtOH 15 vol. %) | 50 | -25.52 | 0.07 | -25.42 | 0.10 |
| | Raw sample | 50 | -25.60 | 0.02 | | 0.18 |

Conclusions

A method for Static Headspace Sampling GC-IRMS was optimized and applied for carbon isotope analysis of methanol and ethanol in aqueous matrix. The isotopic data obtained by SHS GC-IRMS analysis of the alcohols standard mixes demonstrate excellent precision and accuracy, including high sensitivity and low detection limits.

The lifespan of the chromatographic column is extended since only the volatile fraction is injected. During this study, a total of 851 injections (including alcohol standards and real samples) were performed on the TR-WAX chromatographic column. By the end of the study, the column showed no signs of damage, no degradation in chromatographic separation performance, no background issues, and there was no evidence of bleeding. Preventively, 5 cm of the column head was cut three times, and no other maintenance or conditioning was necessary.

The SHS GC-IRMS technique eliminates both the need for time-consuming sample preparation, and the use of expensive specific distillation devices. The newly developed methodology substantially reduces the total sample preparation time from more than 5 hours to few minutes. This results in analytical setup simplification and higher automation, making routine SHS GC-IRMS analysis of ethanol in wines and spirits feasible.

References

1. Yamada K, Yoshida N, Calderone G, et al. (2007) Determination of hydrogen, carbon and oxygen isotope ratios of ethanol in aqueous solution at millimole levels, *Rapid Commun Mass Spectrom*, 21:1431–1437.
2. Cabañero A, Recio JL, Rupérez M (2008) Isotope ratio mass spectrometry coupled to liquid and gas chromatography for wine ethanol characterization, *Rapid Commun Mass Spectrom*, 22:3111–3118.

 Learn more at thermofisher.com/gc-isolink

General Laboratory Equipment. Not for diagnostic procedures. © 2025 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. 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 countries. Please consult your local sales representative for details. **TN-003702-EM-EN 0625**