

EPA TO-15 Analysis Using Hydrogen Carrier Gas and the Agilent HydroInert Source

Ambient air testing using cryogen-free thermal desorption and gas chromatography coupled to a single quadrupole mass spectrometer (GC/MS) with hydrogen gas

Authors

Laura Miles, Hannah Calder, and Helen Martin Markes International Ltd. Llantrisant, UK

Tarun Anumol, Amanda McQuay, and Angela Smith Henry Agilent Technologies, Inc. Wilmington, DE, USA

Abstract

Ambient air monitoring is commonly analyzed with gas chromatography/mass spectrometry (GC/MS) and helium carrier gas. Recent pressure on the helium supply has required organizations to actively investigate hydrogen carrier gas, but most GC/MS analyses have reduced sensitivity and hydrogenation or dechlorination in the sources. This application note describes the use of hydrogen carrier gas and the Agilent Hydrolnert source for GC/MS analysis of humidified canister "air toxics" samples at 100% relative humidity (RH), using cryogen-free systems for thermal desorption preconcentration. Detection of 65 target compounds ranging in volatility from propene to naphthalene is demonstrated, with excellent peak shape and performance well within the criteria set out in U.S. Environmental Protection Agency (EPA) method Toxic Organics-15 (TO-15), including method detection limits (MDLs) as low as 11 parts per trillion by volume (pptv).

Introduction

Monitoring of chemicals in ambient air is necessary to determine the effect they have on the environment and global climate. This monitoring has driven the development of several national and international regulations, primarily in response to increased concern over potentially hazardous volatile organic compounds (VOCs) in ambient (primarily urban) air, industrial emissions, and landfill gas.

Analysis of these VOCs is carried out in accordance with a few standard methods, which require the use of either sorbent tubes (pumped or passive), canisters, or online techniques. Each method has its own advantages and range of applicability, with canister sampling being most popular in the U.S. and China. To achieve the required detection limits using this approach, preconcentration is required to focus analytes and to selectively eliminate bulk constituents. This approach is mandated within the most popular standard method for canisters, U.S. EPA method TO-15.1

Despite the popularity of canister sampling, traditional canister preconcentration technologies are challenged by the ever-greater range of analytes and concentrations of interest. The range of temperatures and humidities at sampling locations is also an issue. High levels of humidity are difficult because the ingress of water to the analytical instrument can negatively impact analyte response and repeatability, as well as reduce the lifetime of the column and detector

The availability of helium has been a concern for several years, and interest in transitioning to alternative carrier gases such as hydrogen has significantly increased. However, existing MS systems have issues with dechlorination of heavily chlorinated compounds. These issues would alter the mass spectra of a peak in the total ion chromatogram (TIC) and lead to potential misidentification of compounds. A newly designed extractor source called the Agilent HydroInert source for the Agilent 5977B Inert Plus GC/MSD addresses these hydrogen-related issues and helps improve performance with hydrogen carrier gas in GC/MS. The HydroInert source with hydrogen carrier gas retains mass spectral fidelity and allows users to continue using existing heliumbased mass spectral libraries and quantitative methods.

This application note shows how the use of a canister autosampler, an innovative trap-based water removal device, and thermal desorption-gas chromatography/mass spectrometry (TD) GC/MS together with hydrogen carrier gas and the Hydrolnert source, allow the analysis of a range of volatile "air toxics" from canisters at 100% RH, in accordance with U.S. EPA method TO-15. Note that, although the term "TO-15" is used by some to describe canister sampling in general, the focus of this work will be on compliance with the specific requirements of the method.

Overview of U.S. EPA method TO-15

The key operations are summarized below.

1. **Sampling:** After cleaning and evacuating the canister, it is brought to the sampling site. The canister valve is opened, and a flow controller draws air through a filter into the canister. After the sampling time corresponding to the set constant flow rate is reached, the canister valve is closed and sealed with caps.

- 2. **Storage:** The sample is kept at ambient temperature and should be analyzed as soon as possible and no later than 20 days after sampling.
- 3. Sample analysis: A known volume of sample is directed from the canister, which is connected to the canister autosampler through a water removal unit and into the multisorbent focusing trap within the concentrator system. The water removal unit will remove most of the water from the sample, and any water vapor remaining in the sample can further be reduced through purging the trap. After the concentration and drying steps are completed, the VOCs are thermally desorbed, entrained in a carrier gas stream, and transferred onto a GC column for separation.
- 4. Compound identification and **quantitation:** Method TO-15 uses GC/MS for qualitative and quantitative analysis of samples. For linear quadrupole MS, monitoring of a wide m/z range (scan mode), or ion-selective scanning (SIM mode) patterns can be used to monitor the relevant target compounds. The mass spectra of the individual peaks in the TIC are examined, and VOCs are identified based on the intensities of quantifier and qualifier ions. The acquired mass spectra are then compared with library spectra (taken under similar conditions) to identify the compound. For any given compound, the abundance of the quantifier ion is compared to its abundance for the compound at known concentrations to determine the concentration of the compound in the sample.

Experimental

Instrumentation

The analytical system used for this study was a multigas CIA Advantage-xr canister autosampler with a Kori-xr water removal device and UNITY-xr thermal desorber, coupled to an Agilent 8890B GC and 5977B single quadrupole GC/MSD system with a HydroInert EI source and 6 mm lens (part number G3870-20448).

Tables 1 and 2 display the canister, TD, GC, and MS parameters.

Standard preparation

A 1 part per million (ppm) standard containing 65 "air toxics" compounds at 1 ppm was diluted in 6 L canisters with nitrogen balance gas to 10 parts per billion by volume (ppbv), unless otherwise stated. RH of 100% was achieved by injecting an appropriate volume of water into the canister.

Results and discussion

A listing of results is provided in Table A1 (see Appendix).

Chromatography

Figure 1 shows typical analyses of the 10 ppbv TO-15 standard at 100% RH, and Figure 2 shows extracted ion chromatograms (EICs) for 16 components spanning the volatility range. Note the excellent peak shape, especially for lighter VOCs, which demonstrates the effectiveness of the Kori-xr module at removing water before analyte trapping at 100% RH.

Another noteworthy aspect is that hydrogen carrier gas decreases the GC cycle time by 40% (45 to 27 minutes), as indicated by the shift in naphthalene from 38 to 23 minutes, while maintaining retention time order for all compounds. Faster chromatography can significantly increase sample throughput.

Table 1. GC and TD parameters.

| Parameter | Value | | | | |
|------------------------------|--|--|--|--|--|
| Gas Chromatograph | Agilent 8890B GC | | | | |
| Column | Agilent J&W DB-624, 60 m × 0.25 mm, 1.40 μm (p/n 123-1364) | | | | |
| Inlet | Splitless | | | | |
| Inlet Temperature | 120 °C | | | | |
| Oven Temperature Program | 30 °C (3 min) 8.3 °C/min to 230 °C (0 min) | | | | |
| Total Run Time | 27 min | | | | |
| MS Transfer Line Temperature | 230 °C | | | | |
| Injection Volume | NA | | | | |
| Carrier Gas | Hydrogen, 2.0 mL/min constant flow | | | | |
| Canister Sampling | | | | | |
| Instrument | CIA Advantage-xr (Markes International) | | | | |
| Sample Volume | Up to 400 mL (for samples of 50 to 100% RH) | | | | |
| | Water Removal | | | | |
| Instrument | Kori-xr (Markes International) | | | | |
| Trap Temperatures | -30 °C/+300 °C | | | | |
| | TD | | | | |
| Instrument | UNITY-xr (Markes International) | | | | |
| Flow Path | 120 °C | | | | |
| Standby Split | 10 mL/min | | | | |
| Sample Flow | 50 mL/min | | | | |
| Trap Purge | 1.0 min at 50 mL/min | | | | |
| Trap Desorption | 2.0 min at 4 mL/min split flow | | | | |
| Cold Trap | Focusing trap: Air Toxics Analyzer (p/n U-T15ATA-2S) | | | | |

Table 2. MS parameters.

| Parameter | Value | | |
|------------------------|----------------------------|--|--|
| Source | HydroInert source | | |
| Mode | Electron ionization, 70 eV | | |
| Source Temperature | 300 °C | | |
| Quadrupole Temperature | 200 °C | | |
| Scan Range | m/z 30 to 300 | | |

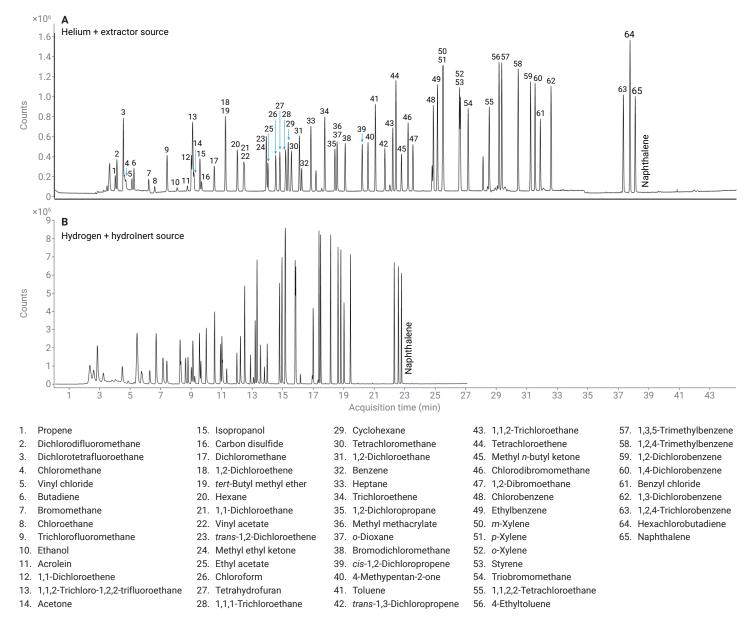


Figure 1. Analysis of 400 mL of a 10 ppbv 65-component TO-15 standard at 100% RH using helium (A) and hydrogen (B) gas. Naphthalene's retention time shifts from 38 to 23 minutes, indicating a 40% reduction in GC cycle time.

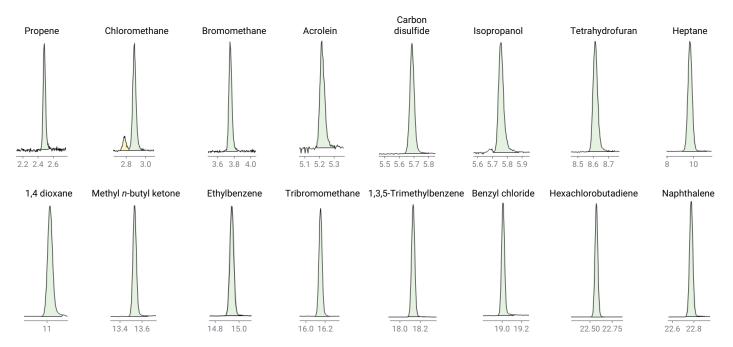


Figure 2. Excellent peak shape of EICs for 16 selected compounds at 2.5 ppbv analyzed using hydrogen gas.

Spectral fidelity

The HydroInert source maintains spectral fidelity by preventing hydrogenation and other reactions that can occur using hydrogen carrier gas. Library match scores (LMS) for all analytes in the 65-component mix were well above 90%, indicating that unwanted source reactions were prevented. Figure 3 shows two examples of high match scores to the National Institute of Standards and Technology helium library (NIST20).

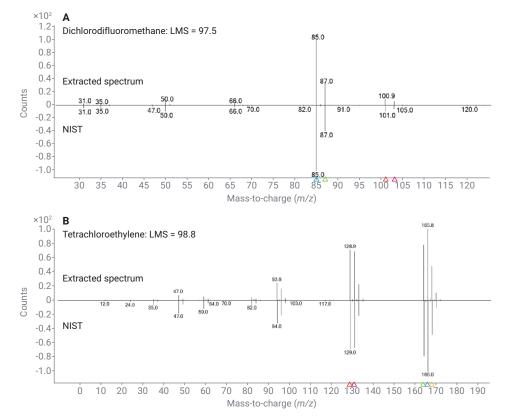


Figure 3. LMS and comparison of extracted versus NIST spectra for dichlorodifluoromethane (A) and tetrachloroethylene (B). Spectral fidelity is maintained.

Linearity

Linearities concerning concentration were calculated at 0.5, 1.25, 2.5, 5, 7.5, and 10 ppbv.

Excellent system linearities were obtained at 100% RH (Table A1), with a mean R² value of 0.999 from 0.50 to 10 ppbv. Figure 4 shows a linearity plot for the 100% RH sample, for the set of 14 compounds spanning the volatility range.

Method detection limits

MDLs were calculated based on seven replicate samples at 0.1 ppbv. 2 To comply with Method TO-15, MDLs are required to be \leq 0.5 ppbv.

Calculations using hydrogen carrier gas gave a mean MDL of 28 pptv for the 28 compounds analyzed (Table 3), which is well within method criteria confirming that method compliance can be achieved, and exceeded, for TO-15 using the HydroInert source. Values ranged from 11 pptv for 4-ethyltoluene, to 53 pptv for carbon disulfide, with a solitary outlier at 113 pptv for propene. These values are all much lower than the requirement of ≤0.5 ppbv.

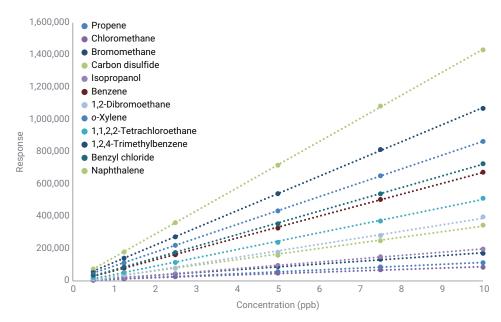


Figure 4. Linearities from 0.5 to 10 ppbv for 14 compounds spanning the range of volatilities, from the 100% RH sample.

Table 3. MDL values for 28 compounds at 0.1 ppbv from the 100% RH sample.

| No. | Compound | Hydrogen MDL (pptv) | |
|-----|---|------------------------|--|
| 1 | Propene | 113 | |
| 2 | Dichlorodifluoromethane | 38 | |
| 3 | Vinyl chloride | 29 | |
| 4 | Butadiene | 33 | |
| 5 | 1,1-Dichloroethene | 24 | |
| 6 | 1,1,2-Trichloro-1,2,2- trifluoroethane | 25 | |
| 7 | Carbon disulfide | 53 | |
| 8 | Dichloromethane | 29 | |
| 9 | tert-Butyl methyl ether | 16 | |
| 10 | Vinyl acetate | 23 | |
| 11 | Methyl ethyl ketone | 42 | |
| 12 | Tetrachloromethane | 17 | |
| 13 | Benzene | 19 | |
| 14 | Heptane | 13 | |

| No. | Compound | Hydrogen MDL (pptv) |
|-----|---------------------------|------------------------|
| 15 | Methyl methacrylate | 34 |
| 16 | Bromodichloromethane | 34 |
| 17 | Toluene | 14 |
| 18 | Tetrachloroethene | 13 |
| 19 | 1,2-Dibromoethane | 28 |
| 20 | Chlorobenzene | 14 |
| 21 | Ethylbenzene | 14 |
| 22 | Styrene | 23 |
| 23 | 1,1,2,2-Tetrachloroethane | 16 |
| 24 | 4-Ethyltoluene | 11 |
| 25 | 1,3,5-Trimethylbenzene | 14 |
| 26 | 1,2-Dichlorobenzene | 17 |
| 27 | Hexachlorobutadiene | 23 |
| 28 | Naphthalene | 50 |
| | Average | 28 |

Reproducibility

Method TO-15 requires that the calculated relative standard deviations (RSDs) for the relative response factors (RRFs) for each compound in the calibration table must be less than 30%, with at most two exceptions up to a

limit of 40%. Results were well within the requirements of method TO-15 for 100% RH at 7.47% RSD. Further, for 10 replicates of 10 ppbv at 100% RH, average area RSD was 1.22%. For 50 injections of varying concentrations of 100% RH, average retention time

RSD was 0.09%, where criteria state <1% variation (Figure 5). Efficient water management using UNITY-Kori-CIA Advantage-xr enables stable retention times and highly reproducible peak area responses, even with hydrogen carrier gas.

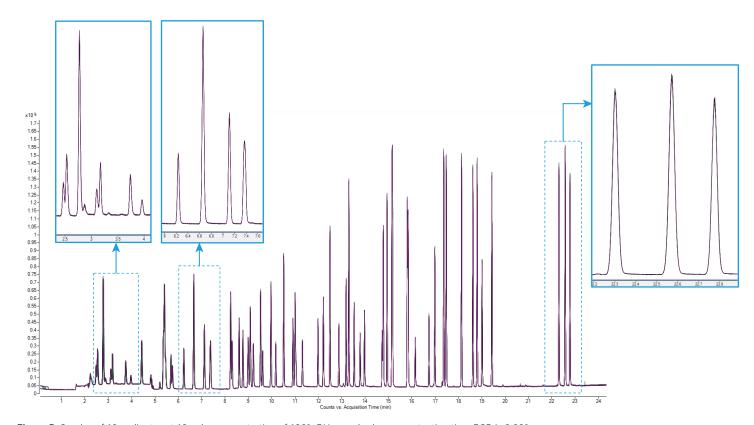
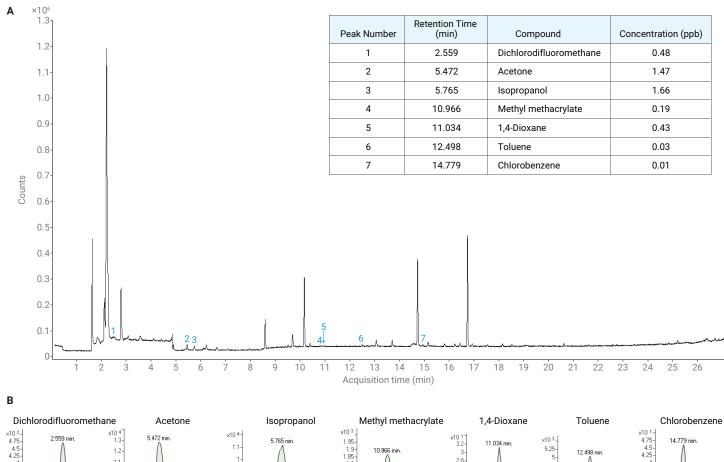


Figure 5. Overlay of 10 replicates at 10 ppbv concentration of 100% RH sample. Average retention time RSD is 0.09%.

Real air sample

To illustrate the performance of the system for a real air sample, 400 mL of lab air was analyzed under the same conditions as described previously. Seven components from the 65-member TO-15 list were found to be at quantifiable levels (Figure 6).



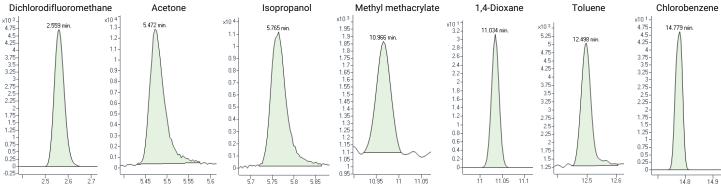


Figure 6. (A) Real sample chromatogram generated from analysis of 400 mL of rural air, using the conditions previously described. Compounds from the TO-15 listing are indicated. (B) Zoom of seven compounds detected from the TO-15 list.

Conclusion

The Agilent 8890 GC coupled to the Agilent 5977B single quadrupole MS and CIA Advantage—Kori-xr—UNITY-xr preconcentration system with hydrogen carrier gas allows confident analysis of "air toxics" in humid environments, in accordance with U.S. EPA method TO-15.

Key features of the results are the excellent chromatographic performance for the analysis of a 65-component T0-15 mix (from propene to naphthalene) at 100% relative humidity. Performance was well within the requirements of method T0-15, with method detection limits as low as 11 pptv.

Crucially, this performance has been achieved for even the most volatile of components in the TO-15 mix, due to the efficient and selective removal of water from humid air streams without compromising the analysis of VOCs or polar species. In addition, the system uses electrical trap cooling, eliminating the cost and inconvenience associated with liquid cryogen. The Agilent HydroInert source with hydrogen carrier gas retained mass spectral fidelity, allowing users to continue the use of existing helium-based mass spectral libraries and quantitative methods. The robustness and reliability of the GC/MSD allows long-term operation of the system while generating data in compliance with U.S. EPA TO-15 requirements.

References

- Compendium Method TO-15:
 Determination of Volatile Organic Compounds (VOCs) in Air Collected in Specially-Prepared Canisters and Analyzed by Gas Chromatography/ Mass Spectrometry (GC/MS), Compendium of Methods for the Determination of Toxic Organic Compounds in Ambient Air (second edition), US EPA 1999.
 https://19january2017snapshot.epa.gov/homeland-security-research/epa-air-method-toxic-organics-15-15-determination-volatile-organic_.html
- Ambient Air Determination of Volatile Organic Compounds Tank Sampling/Gas Chromatography-Mass Spectrometry, Ministry of Ecology and Environment of the People's Republic of China (mee. gov.cn). https://www.mee.gov.cn/ ywgz/fgbz/bz/bzwb/jcffbz/201510/ t20151030_315940.shtml

Appendix

Many conventional systems for canister analysis use liquid cryogen to trap VOCs. Markes' systems instead use electrical (Peltier) cooling, allowing quantitative retention of the most VOCs from large sample volumes, without incurring the cost of liquid cryogen.

Method TO-15 states that any canister that has not tested clean (compared to direct analysis of humidified zero air of less than 0.2 ppbv of targeted VOCs) should not be used.

MDLs were calculated based on 99% confidence for seven values (MDL = $3.143 \times \text{standard deviation} \times \text{concentration}$).

Table A1. Data comparison of helium and hydrogen carrier gas results obtained for TO-15 standards at 100% RH. Linearity (R^2) values were generated for the concentration range 0.22 to 10 ppbv for analysis with helium carrier gas, and 0.5 to 10 ppv for hydrogen carrier gas.

| | | Helium Carrier (100% RH) | | | Hydrogen Carrier (100% RH) | | | |
|-----|---------------------------------------|--------------------------|-------------|-------------|----------------------------|-------------|-------------|--|
| No. | Compound | RT (min) | RRF RSD (%) | RRF RSD (%) | RT (min) | RRF RSD (%) | RRF RSD (%) | |
| 1 | Propene | 4.894 | 0.9997 | 6.3 | 2.47 | 0.9994 | 12.4% | |
| 2 | Dichlorodifluoromethane | 5.032 | 0.9998 | 5.6 | 2.53 | 1.0000 | 8.4% | |
| 3 | Dichlorotetrafluoroethane | 5.500 | 0.9997 | 7.7 | 2.77 | 0.9999 | 4.2% | |
| 4 | Chloromethane | 5.686 | 0.9808 | 11.0 | 2.88 | 0.9929 | 9.3% | |
| 5 | Vinyl chloride | 6.122 | 0.9994 | 4.6 | 3.10 | 0.9997 | 10.1% | |
| 6 | Butadiene | 6.276 | 0.9998 | 2.0 | 3.17 | 0.9999 | 12.4% | |
| 7 | Bromomethane | 7.346 | 0.9988 | 14.4 | 3.74 | 0.9988 | 10.3% | |
| 8 | Chloroethane | 7.723 | 0.9870 | 9.1 | 3.95 | 0.9999 | 5.2% | |
| 9 | Trichlorofluoromethane | 8.646 | 0.9999 | 6.2 | 4.42 | 0.9996 | 5.6% | |
| 10 | Ethanol | 9.299 | 0.9997 | 24.4 | 4.80 | 0.9990 | 20.9% | |
| 11 | Acrolein | 9.925 | 0.9993 | 9.3 | 5.20 | 0.9988 | 18.9% | |
| 12 | 1,1-Dichloroethene | 10.258 | 0.9998 | 1.2 | 5.35 | 0.9999 | 13.2% | |
| 13 | 1,1,2-Trichloro-1,2,2-trifluoroethane | 10.337 | 1.0000 | 4.8 | 5.40 | 0.9998 | 2.2% | |
| 14 | Acetone | 10.401 | 1.0000 | 2.7 | 5.45 | 0.9999 | 5.1% | |
| 15 | Isopropanol | 10.868 | 0.9981 | 18.8 | 5.74 | 0.9996 | 14.2% | |
| 16 | Carbon disulfide | 10.884 | 0.9999 | 0.9 | 5.67 | 0.9984 | 5.6% | |
| 17 | Dichloromethane | 11.657 | 0.9998 | 2.0 | 6.23 | 0.9995 | 8.7% | |
| 18 | 1,2-Dichloroethene | 12.461 | 0.9999 | 1.6 | 6.65 | 0.9997 | 11.8% | |
| 19 | tert-Butyl methyl ether | 12.513 | 0.9997 | 4.2 | 6.65 | 0.9999 | 3.4% | |
| 20 | Hexane | 13.285 | 0.9956 | 13.6 | 7.11 | 0.9999 | 2.3% | |
| 21 | 1,1-Dichloroethane | 13.578 | 1.0000 | 5.5 | 7.36 | 0.9996 | 10.1% | |
| 22 | Vinyl acetate | 13.737 | 0.9998 | 1.5 | 7.39 | 0.9981 | 12.5% | |
| 23 | trans-1,2-Dichloroethene | 15.112 | 0.9998 | 3.6 | 8.24 | 0.9997 | 11.4% | |
| 24 | Methyl ethyl ketone | 15.127 | 0.9998 | 9.2 | 8.23 | 0.9998 | 6.9% | |
| 25 | Ethyl acetate | 15.314 | 0.9999 | 5.5 | 8.30 | 1.0000 | 4.7% | |
| 26 | Chloroform | 15.904 | 0.9999 | 5.5 | 8.76 | 0.9984 | 8.6% | |
| 27 | Tetrahydrofuran | 15.912 | 0.9998 | 9.2 | 8.60 | 0.9997 | 6.8% | |
| 28 | 1,1,1-Trichloroethane | 16.447 | 0.9999 | 8.2 | 8.99 | 0.9969 | 7.9% | |
| 29 | Cyclohexane | 16.637 | 0.9999 | 8.8 | 9.08 | 0.9997 | 2.6% | |
| 30 | Tetrachloromethane | 16.902 | 1.0000 | 7.4 | 9.21 | 0.9942 | 11.1% | |
| 31 | 1,2-Dichloroethane | 17.378 | 1.0000 | 3.5 | 9.61 | 0.9999 | 5.6% | |
| 32 | Benzene | 17.390 | 0.9999 | 6.1 | 9.52 | 0.9998 | 2.0% | |
| 33 | Heptane | 18.075 | 0.9995 | 19.0 | 9.97 | 0.9998 | 2.8% | |
| 34 | Trichloroethene | 19.022 | 0.9999 | 5.1 | 10.51 | 0.9990 | 10.0% | |
| 35 | 1,2-Dichloropropane | 19.557 | 0.9999 | 7.1 | 10.91 | 0.9990 | 4.9% | |
| 36 | Methyl methacrylate | 19.822 | 0.9989 | 2.9 | 11.00 | 0.9997 | 2.8% | |
| 37 | <i>p</i> -Dioxane | 19.914 | 0.9998 | 12.6 | 11.01 | 0.9981 | 15.6% | |
| 38 | Bromodichloromethane | 20.227 | 0.9999 | 6.2 | 11.32 | 0.9953 | 11.9% | |
| 39 | cis-1,3-Dichloropropene | 21.399 | 1.0000 | 4.5 | 11.98 | 0.9961 | 12.9% | |
| 40 | 4-Methylpentan-2-one | 21.760 | 0.9999 | 2.8 | 12.21 | 0.9996 | 5.6% | |
| 41 | Toluene | 22.326 | 0.9999 | 16.3 | 12.50 | 1.0000 | 4.1% | |
| 42 | trans-1,3-Dichloropropene | 22.810 | 0.9997 | 2.9 | 12.88 | 0.9973 | 13.1% | |
| 43 | 1,1,2-Trichloroethane | 23.305 | 1.0000 | 7.9 | 13.19 | 0.9990 | 6.3% | |

| | | Helium Carrier (100% RH) | | Hydrogen Carrier (100% RH) | | | |
|------------|-------------------------------|--------------------------|-------------|----------------------------|----------|-------------|-------------|
| No. | Compound | RT (min) | RRF RSD (%) | RRF RSD (%) | RT (min) | RRF RSD (%) | RRF RSD (%) |
| 44 | Tetrachloroethene | 23.828 | 1.0000 | 8.8 | 13.31 | 0.9996 | 2.9% |
| 45 | Methyl <i>n</i> -butyl ketone | 23.959 | 0.9998 | 2.5 | 13.53 | 0.9942 | 17.0% |
| 46 | Chlorodibromomethane | 24.398 | 0.9999 | 4.7 | 13.79 | 0.9973 | 12.6% |
| 47 | 1,2-Dibromoethane | 24.735 | 1.0000 | 4.4 | 13.98 | 0.9973 | 14.2% |
| 48 | Chlorobenzene | 26.102 | 1.0000 | 11.9 | 14.78 | 0.9998 | 2.2% |
| 49 | Ethylbenzene | 26.407 | 0.9999 | 20.2 | 14.94 | 1.0000 | 4.6% |
| 50 | m-Xylene | 26.732 | 1.0000 | 20.8 | 15.16 | 0.9999 | 4.7% |
| 51 | <i>p</i> -Xylene | 26.732 | 1.0000 | 20.8 | 15.16 | 0.9999 | 4.7% |
| 52 | o-Xylene | 27.837 | 1.0000 | 25.2 | 15.81 | 1.0000 | 5.7% |
| 53 | Styrene | 27.857 | 0.9999 | 10.3 | 15.84 | 1.0000 | 4.1% |
| 54 | Tribromomethane | 28.376 | 0.9998 | 4.4 | 16.15 | 0.9976 | 6.8% |
| 5 5 | 1,1,2,2-Tetrachloroethane | 29.624 | 0.9999 | 6.8 | 16.99 | 0.9990 | 8.1% |
| 56 | 4-Ethyltoluene | 30.385 | 0.9999 | 6.3 | 17.37 | 1.0000 | 4.1% |
| 57 | 1,3,5-Trimethylbenzene | 30.551 | 1.0000 | 19.6 | 18.13 | 1.0000 | 4.5% |
| 58 | 1,2,4-Trimethylbenzene | 31.653 | 1.0000 | 10.4 | 17.48 | 0.9999 | 4.6% |
| 59 | 1,2-Dichlorobenzene | 32.485 | 0.9999 | 3.9 | 18.62 | 1.0000 | 1.0% |
| 60 | 1,4-Dichlorobenzene | 32.738 | 0.9999 | 3.3 | 18.80 | 0.9999 | 2.4% |
| 61 | Benzyl chloride | 33.107 | 0.9998 | 2.3 | 19.01 | 0.9998 | 2.0% |
| 62 | 1,3-Dichlorobenzene | 33.840 | 0.9999 | 7.4 | 19.43 | 0.9999 | 3.8% |
| 63 | 1,2,4-Trichlorobenzene | 38.594 | 0.9965 | 18.9 | 22.30 | 0.9998 | 2.2% |
| 64 | Hexachlorobutadiene | 39.121 | 0.9997 | 9.4 | 22.57 | 0.9988 | 7.5% |
| 65 | Naphthalene | 39.315 | 0.9975 | 19.7 | 22.78 | 1.0000 | 1.6% |
| | Mean values | | 0.9992 | 8.5 | | 0.9990 | 7.5% |

www.agilent.com

DE14855333

This information is subject to change without notice.

