



Application Note 059

Direct Desorption of Car Trim Materials for VOC and SVOC Analysis in Accordance with VDA Method 278

Application Note



MARKES
international

Abstract

In this application note, we demonstrate the suitability of Markes' TD-100 automated thermal desorption system for performing the analysis of volatile and semivolatile organic compounds (VOCs and SVOCs) in three types of automotive trim materials (polypropylene, artificial leather, and foam) in accordance with the German Association of the Automotive Industry (VDA) Method 278.

Introduction

The German Association of the Automotive Industry (VDA) consists of automobile manufacturers, their development partners, and their suppliers. The VDA promotes German motor trade interest by participating in the development of economic, transport, and environmental policy, and setting appropriate technical standards. Given the international nature of the automotive industry, VDA actions have an impact on subsidiaries and suppliers based in other countries.

Concern is growing worldwide over the quality of air in vehicle cabins. Many people, including vulnerable groups such as children and the elderly, spend a significant proportion of their time in cars and are at risk of being exposed to odorous or toxic organic chemicals that may be emitted from car interior trim components. The levels of volatile and semivolatile organic compounds (VOCs and SVOCs) found in car cabin air can be several times higher than normal ambient air levels (indoor or outdoor) – see Application note 033.



Agilent Technologies

In response to (S)VOC concerns, the VDA developed a series of methods for assessing (S)VOC emissions from car trim components such as upholstery, carpets, glues, sealing compounds, plastics, foam, leather, coatings, and paint. One such test, known as VDA Method 278, requires the assessment of two values:

- **VOC value:** The sum of the VOCs and SVOCs up to n-C₂₅ emitted under one set of specified analytical conditions.
- **FOG value:** The sum of the SVOCs (from n-C₁₄ to n-C₃₂) emitted under a different set of specified test conditions.

VDA Method 278 specifies direct thermal desorption (TD) (using heat and a flow of inert gas) of small samples of relatively homogeneous materials used in car trim components (leather, plastic, and so forth). VOCs and SVOCs are extracted from the sample into the gas stream and are refocused onto a secondary focusing trap prior to injection into a GC/MS or GC/FID system for analysis.

VDA Method 278 comprises two extraction stages:

- **VOC analysis:** This involves desorbing the sample at 90 °C for 30 minutes to extract compounds with volatilities up to n-C₂₅. This is followed by semiquantitative analysis of each compound as micrograms of toluene equivalents per gram of sample.
- **FOG analysis:** This involves desorbing the sample at 120 °C for 60 minutes to extract compounds ranging in volatility from n-C₁₄ to n-C₃₂. This is followed by semiquantitative analysis of each compound as micrograms of *n*-hexadecane equivalents per gram of sample.

Experimental

Samples

Polypropylene: 30 ±5 mg
 Leather: 10 ±3 mg
 Foam: 15 ±5 mg

All samples were weighed directly into an empty unfritted glass sample tube with restrictions at either 15 mm or 30 mm, depending on the sample material.

A calibration solution of toluene and hexadecane in methanol at a concentration of 500 ng/μL was introduced into the sampling end of a sample tube packed with 200 mg of Tenax TA sorbent in a stream of pure inert gas.

Step 1: Determination of VOC value

Parameter	Value
Instrument:	TD-100 (Markes International)
Desorption temperature:	90 °C
Desorption time:	30 minutes
Focusing trap:	General-purpose hydrophobic (Tenax TA–graphitised carbon black)
Focusing trap low:	–30 °C
Focusing trap high:	300 °C
Focusing trap hold:	3 minutes
Flow path temperature:	200 °C
Trap flow:	20 mL/min
Split flow (inlet and outlet):	20 mL/min
Carrier gas flow:	1.3 mL/min (constant flow)
Temperature program:	40 °C (2 minutes), then 3 °C/min to 92 °C, then 5 °C/min to 160 °C, then 10 °C/min to 280 °C (10 minutes)
Mass range:	<i>m/z</i> 29-280
Auxiliary temperature:	280 °C

Step 2: Determination of FOG value

Parameter	Value
Instrument:	TD-100 (Markes International)
Desorption temperature:	120 °C
Desorption time:	60 minutes
Focusing trap:	General-purpose hydrophobic (Tenax TA–graphitised carbon black)
Focusing trap low:	–30 °C
Focusing trap high:	300 °C
Focusing trap hold:	5 minutes
Flow path temperature:	200 °C
Trap flow:	20 mL/min
Split flow (inlet and outlet):	20 mL/min
Temperature program:	50 °C (2 minutes), then 25 °C/min to 160 °C, then 10 °C/min to 310 °C (30 minutes)
Mass range:	<i>m/z</i> 29-370
Auxiliary temperature:	280 °C

Results and Discussion

VOC and FOG analyses

Results of the VOC and FOG analyses for samples of polypropylene, artificial leather, and foam, using Markes' TD-100 automated thermal desorber, are presented in Figures 1-6.

System inertness and recovery of active compounds

To demonstrate system suitability (inertness) for polar species, an activity test mix containing 18 compounds was prepared at a concentration of $0.4 \mu\text{g}/\mu\text{L} \pm 0.1 \mu\text{g}/\mu\text{L}$ (Table 1).

Table 1. Activity test mix and associated component concentration levels.

No.	Compound	Concentration ($\mu\text{g}/\mu\text{L}$)
1	Benzene	0.44
2	<i>n</i> -Heptane	0.35
3	Toluene	0.43
4	<i>n</i> -Octane	0.30
5	<i>p</i> -Xylene	0.43
6	<i>o</i> -Xylene	0.44
7	<i>n</i> -Nonane	0.36
8	<i>n</i> -Decane	0.37
9	2-Ethylhexan-1-ol	0.42
10	<i>n</i> -Undecane	0.37
11	2,6-Dimethylphenol	0.45
12	<i>n</i> -Dodecane	0.38
13	<i>n</i> -Tridecane	0.38
14	<i>n</i> -Tetradecane	0.38
15	Dicyclohexylamine	0.45
16	<i>n</i> -Pentadecane	0.38
17	<i>n</i> -Hexadecane	0.39
18	Bis(2-ethylhexyl) adipate	0.46

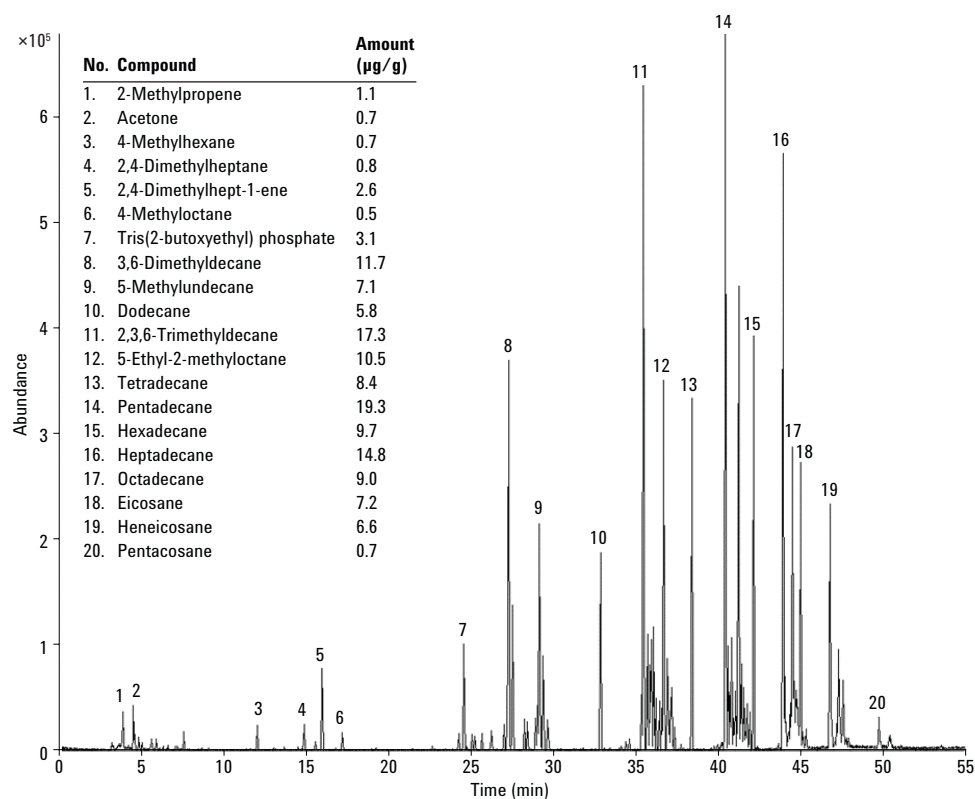


Figure 1. VOC analysis of a 31.4 mg sample of polypropylene.

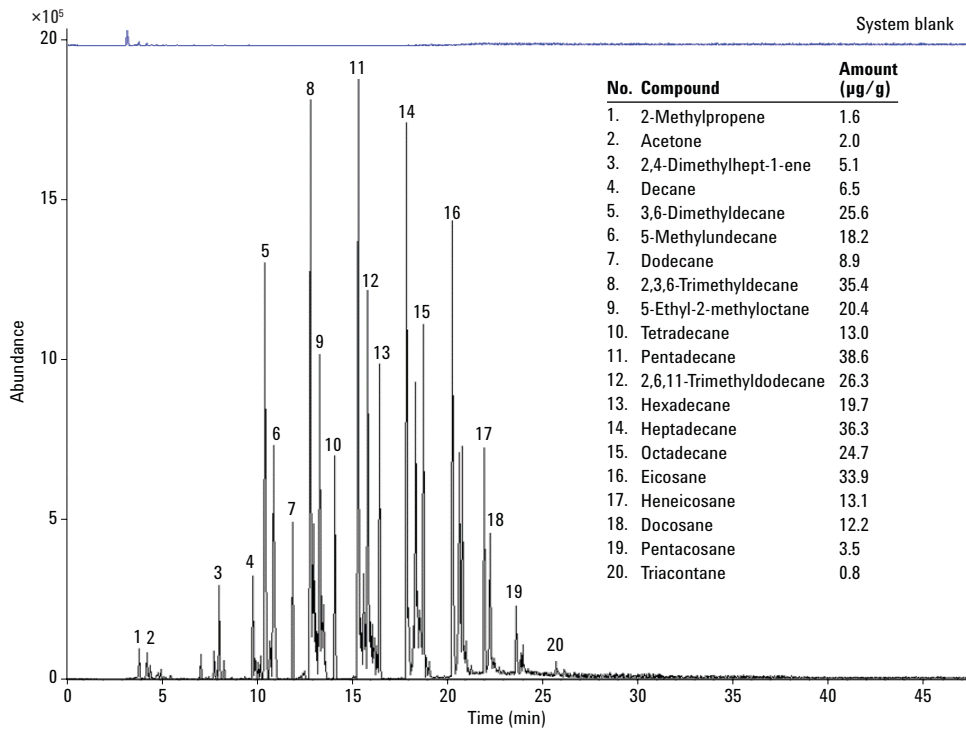


Figure 2. FOG analysis of a sample of polypropylene, followed by a system blank, showing no carryover of high-boiling compounds.

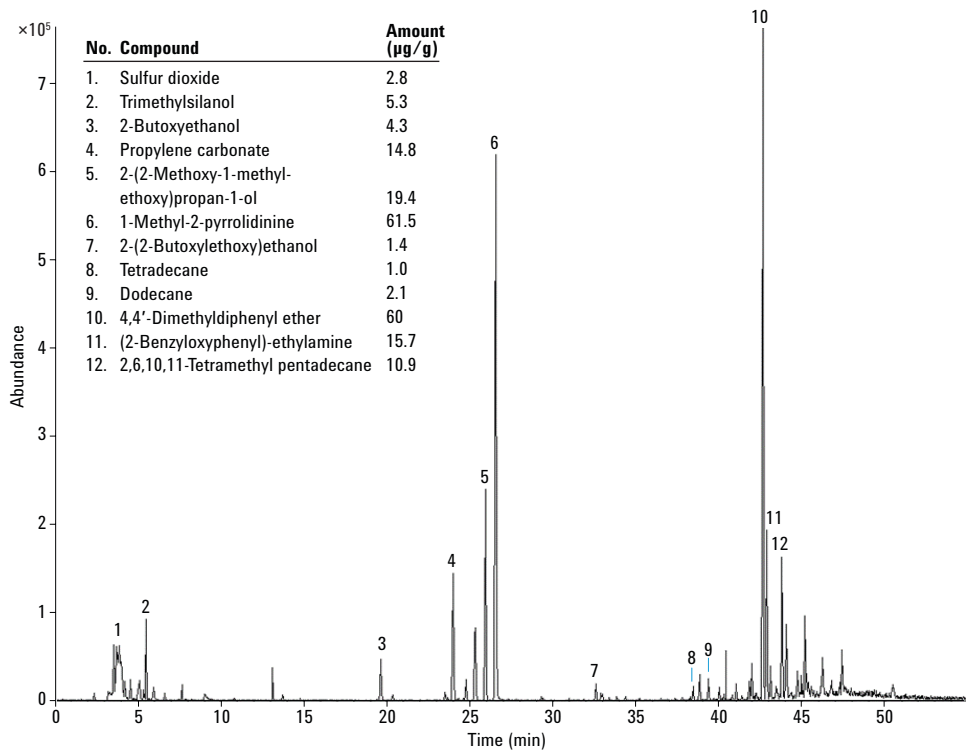


Figure 3. VOC analysis of a 12.5 mg sample of artificial leather.

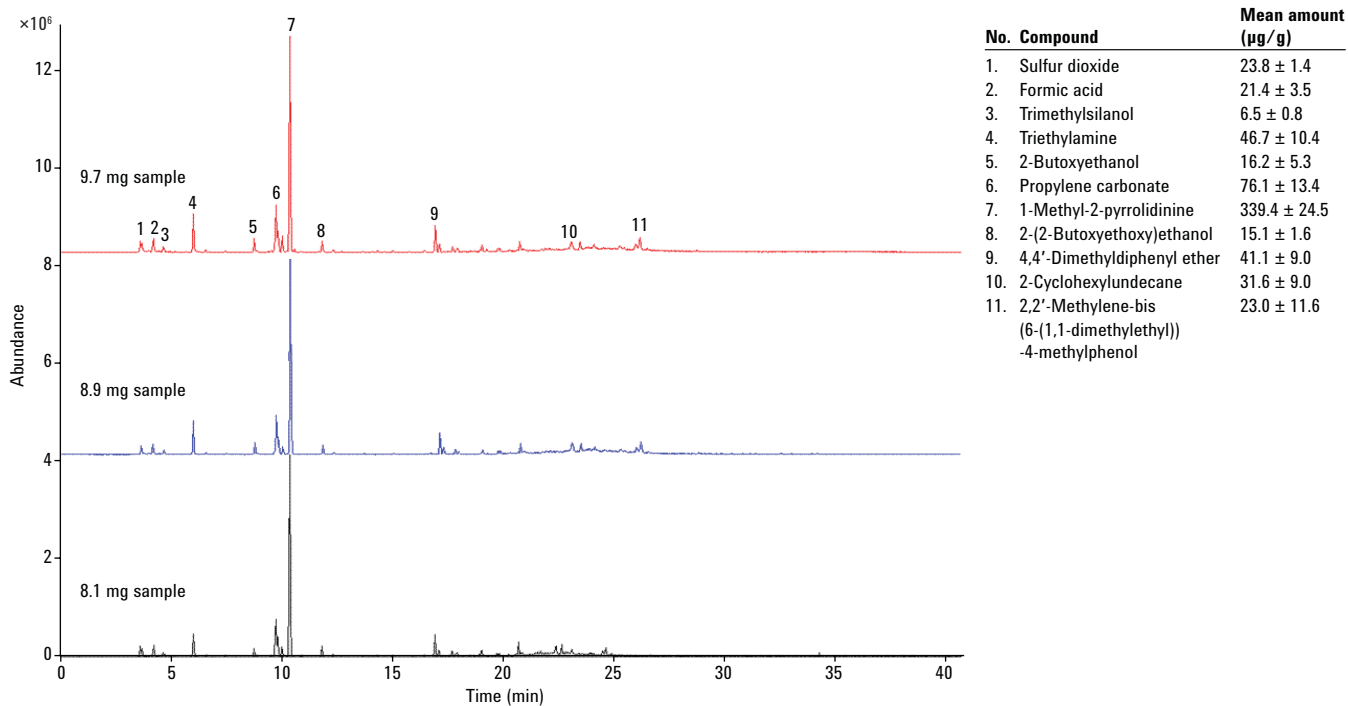


Figure 4. FOG analysis of three samples of artificial leather.

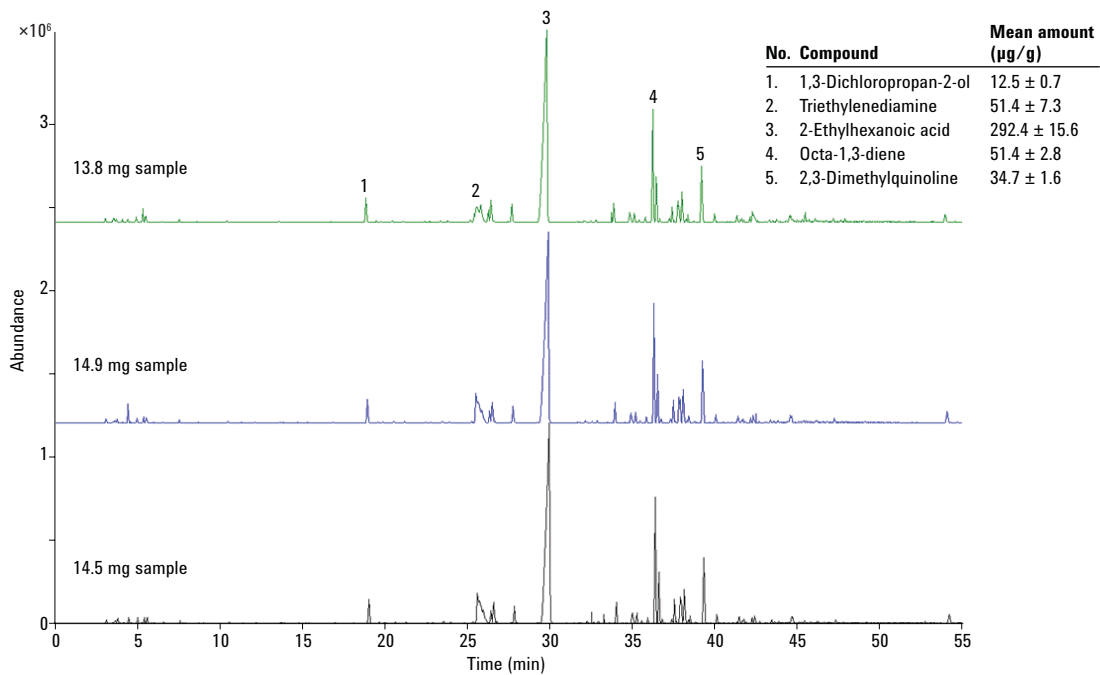


Figure 5. VOC analysis of three samples of foam.

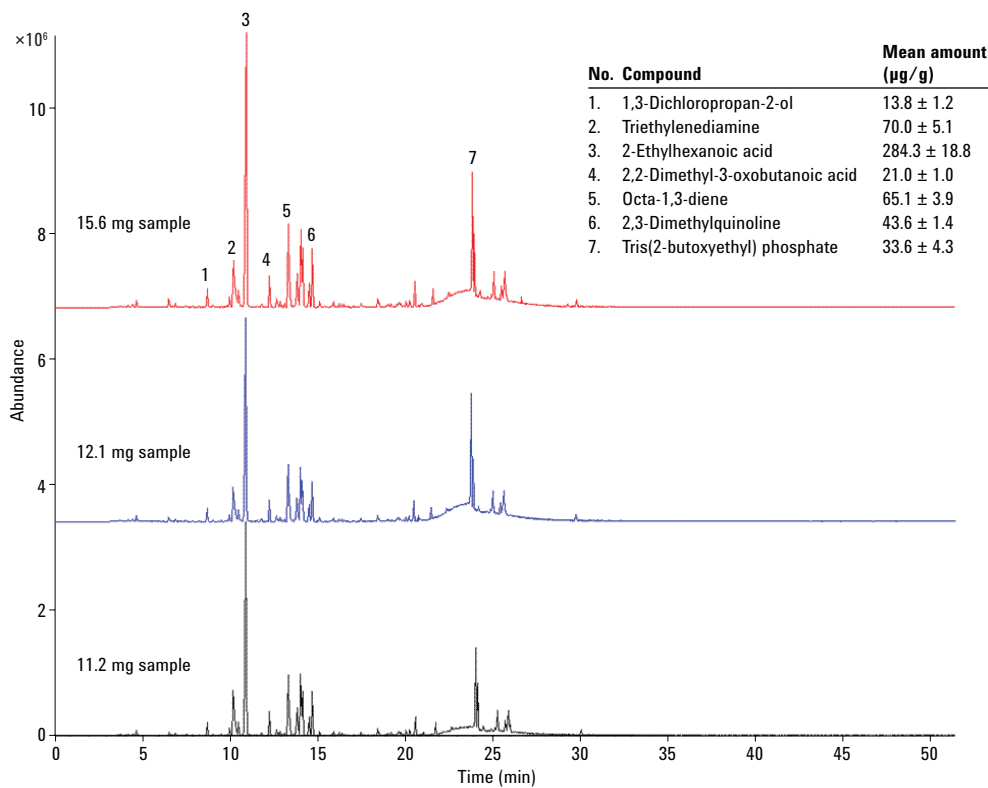


Figure 6. FOG analysis of three samples of foam.

A 1 µL sample of the activity mix was injected into the sampling end of a conditioned Tenax TA tube in a stream of inert gas using the Markes Calibration Solution Loading Rig (CSLR). During analysis, the split portion of the sample was re-collected into a second conditioned Tenax TA tube. Figure 7 shows the chromatogram obtained from the analysis of a Tenax TA tube loaded with 1 µL of activity test mix and desorbed using the VOC analysis conditions. This procedure was repeated a second time, giving three analyses from the one sample (Figure 8).

The lack of bias between the three chromatograms shows that there are no losses in the system, thus validating recovery of all analytes.

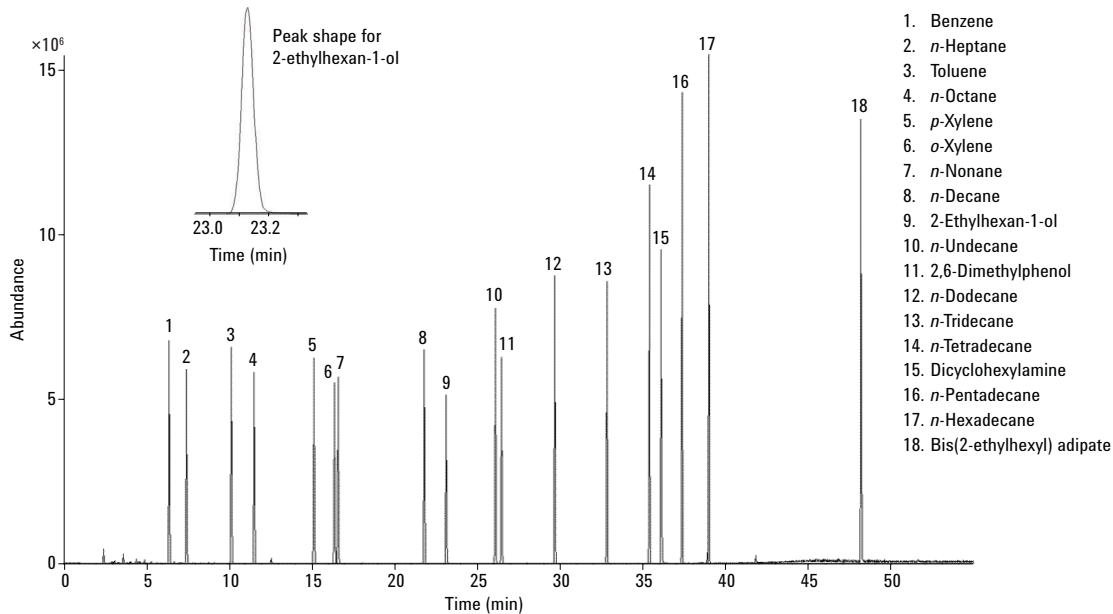


Figure 7. Analysis of 0.4 $\mu\text{g}/\mu\text{L}$ activity test mix spiked onto a Tenax TA tube.

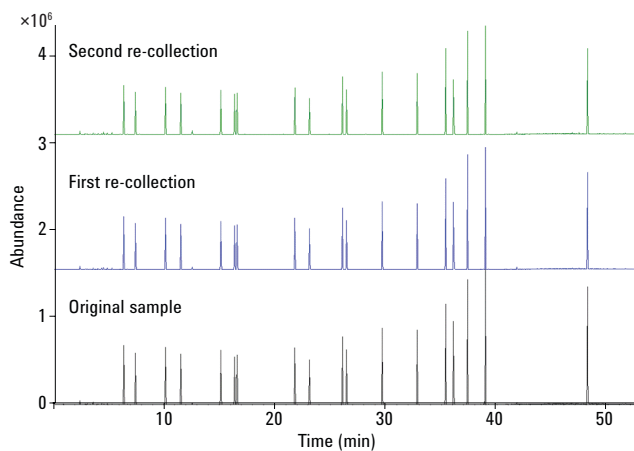


Figure 8. Repeated analyses of the activity test mix.

Repeatability of toluene analysis

A 0.5 μg sample of toluene was injected into 20 tubes packed with Tenax TA. These 20 samples were analyzed for 1 week and produced the following values:

- Average recovery: 100%
- Standard deviation: 5%
- Minimum recovery: 92% (80%)
- Maximum recovery: 110% (120%)

These were found to be within the limits outlined in VDA Method 278 (bracketed).

Linearity and limit of determination/detection

System linearity was demonstrated over the range 0.01-5 μg by preparing a range of liquid standards of toluene and eicosane ($n\text{-C}_{20}$) in methanol, and dotriacontane ($n\text{-C}_{32}$) in *n*-hexane. These were introduced into the sampling end of conditioned Tenax TA tubes in a stream of clean inert gas.

Detection and determination limits were determined in accordance with the procedure described in VDA Method 278. Split and GC conditions used correspond with the VOC and FOG method parameters, respectively. All tubes were desorbed at 320 $^{\circ}\text{C}$ to ensure complete recovery of the standard solutions from the Tenax TA sorbent.

Figures 9-11 show plots of the linearity of the system for toluene, eicosane, and dotriacontane.

In each case, LOD and LOQ values were determined using the 0.01-0.10 μg loading ranges (Table 2).

Table 2. Values for limit of detection (LOD) and limit of determination (LOQ) for the three compounds investigated over the loading range 0.01-0.10 μg .

Compound	LOD*	LOQ
Toluene	0.006 μg (<0.04 μg)	0.017 μg
Eicosane	0.006 μg (<0.06 μg)	0.02 μg
Dotriacontane	0.009 μg (<0.06 μg)	0.02 μg

* The limit cited in VDA Method 278 is given in parentheses.

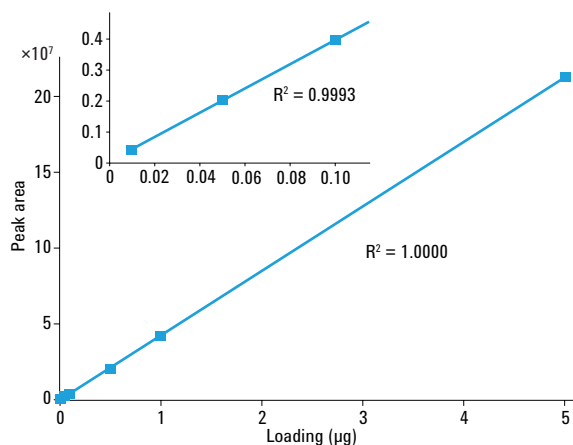


Figure 9. Linearity of toluene over the loading range 0.01-5 μg and (inset) over the range 0.01-0.10 μg .

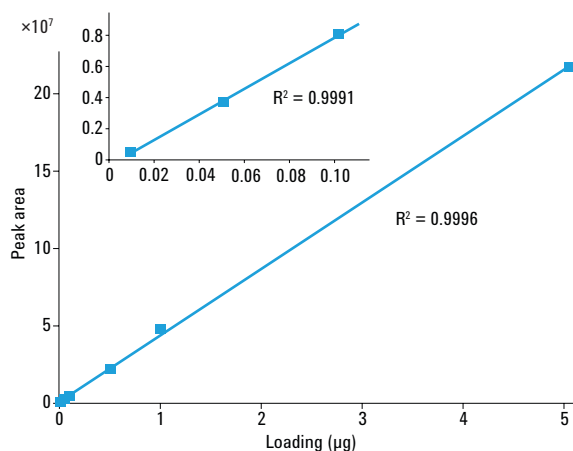


Figure 10. Linearity of eicosane over the loading range 0.01-5 μg and (inset) over the range 0.01-0.10 μg .

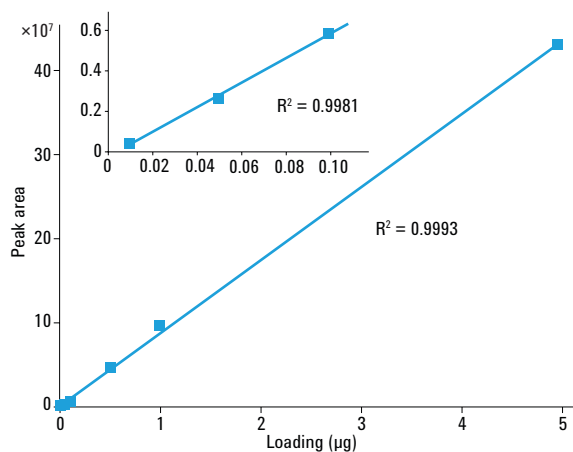


Figure 11. Linearity of dotriacontane over the loading range 0.01-5 μg and (inset) over the range 0.01-0.10 μg .

Conclusions

The suitability of Markes' TD-100 thermal desorber for performing the analysis of automotive trim materials in accordance with VDA Method 278 has been demonstrated.

Analyses of a variety of typical trim materials/standard solutions demonstrated:

- Excellent linearity.
- Low detection/determination levels.
- Excellent peak shape for reactive (polar) compounds, demonstrating good system inertness.
- Repeatable results even for very high-boiling compounds.

Key features of the TD-100 system that are essential for VDA Method 278 include:

- Leak-testing of each tube prior to analysis to ensure data integrity.
- Effective sealing of all sample tubes loaded onto the autosampler before and after analysis to prevent contamination and loss of volatiles.
- Inert flow path for compatibility with reactive species.
- Backflushed and electrically-cooled (cryogen-free) focusing trap for efficient trapping and desorption of compounds from C₂ to n-C₄₄.
- Full EPC control of carrier gas through the TD system for optimum retention time stability and improved qualitative identification of minor compounds.
- Quantitative re-collection for repeat analysis or method validation, to demonstrate quantitative recovery through the system and allow method validation.

Trademarks

CSLR™ and TD-100™ are trademarks of Markes International.

Tenax® is a registered trademark of Buchem B.V..

Applications were performed under the stated analytical conditions. Operation under different conditions, or with incompatible sample matrices, may impact the performance shown.

For More Information

These data represent typical results. For more information on our products and services, visit our Web site at www.agilent.com/chem.

www.agilent.com/chem

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 2017
Printed in the USA
July 12, 2017
5991-8265EN



Agilent Technologies