

# Enhanced Productivity for Residual Solvent Analysis in Pharmaceutical Products According to USP 467 by Using a New Valve-and-Loop Static Headspace Sampler

Xiaoteng Gong<sup>1</sup>, Benjamin Webber<sup>1</sup>, Gayatri Trivedi<sup>1</sup>, Danny Hower<sup>1</sup>, Dujuan Lu<sup>1</sup>, Julian Gulbinski<sup>1</sup>, Cristian I. Cojocariu<sup>2</sup>, Giulia Riccardino<sup>2</sup>, Daniela Cavagnino<sup>3</sup>, Manuela Bergna<sup>3</sup>, Carlos Garcia<sup>4</sup>

<sup>1</sup> SGS, Fairfield, NJ, <sup>2</sup> Thermo Fisher Scientific, Runcorn, UK, <sup>3</sup> Thermo Fisher Scientific, Milan, Italy, <sup>4</sup> Thermo Fisher Scientific, Austin, TX

## ABSTRACT

Residual solvents can be present in pharmaceutical products as a result of the manufacturing process or as a contamination during packaging, warehouse storage or transportation. In order to prevent patients from potentially hazardous effects of those chemicals, pharmaceutical products need to be tested to ensure that any solvent residuals have been efficiently removed during manufacturing processes or, if present, their concentration is compatible with the accepted safety requirements. United States Pharmacopoeia (USP) Method <467> regulates the testing procedure and establishes recommended acceptable levels and the instrument performance criteria. Gas chromatography coupled with valve-and-loop static headspace sampling (GC-HS) is the technique of choice for identification and quantitation of residual solvents in pharmaceutical products. The USP compliance of the new Thermo Scientific™ TriPlus™ 500 Static Headspace Sampler coupled with the Thermo Scientific™ TRACE™ 1310 GC in a routine testing environment is shown here.

## INTRODUCTION

Solvents are widely used in the synthesis of pharmaceutical products, substances and excipients although they cannot always be completely removed during the manufacturing processes. To ensure patients' safety, final products are tested to assess whether the solvents used during the manufacturing processes have been efficiently removed or, if still present, their concentration is within the accepted limits. As organic solvents have relatively low boiling points and are thermally stable the analytical method of choice for Class 1 and Class 2 residual solvent determination is headspace-gas chromatography, with detection using either flame ionization detection (FID) or mass spectrometry (MS). Headspace sampling allows for the extraction of semi-volatile and volatile compounds from complex liquid and solid matrices in a fast and simple way without the need for time-consuming sample preparation. A TriPlus 500 HS autosampler was coupled to a TRACE 1310 GC equipped with a Thermo Scientific™ Instant Connect Split/Splitless SSL Injector and a Thermo Scientific™ Instant Connect FID and used for the determination of residual solvent content in water-soluble and water-insoluble pharmaceuticals according to the United States Pharmacopoeia <467> method (USP).<sup>1</sup>

## MATERIALS AND METHODS

USP <467> Class 1, Class 2A and Class 2B residual solvent solutions in dimethylsulfoxide (DMSO) were purchased from (Restek®, P/N 36279, 36012, 36280 respectively). Stock and standard solutions for procedures A, B, C were diluted in HPLC-MS grade water or GC headspace grade DMSO as reported in the USP <467> method. Dispersive aspirin (acetylsalicylic acid, 75 mg) and common pain relief tablets (paracetamol, 500 mg and caffeine, 65 mg) were used to prepare sample stock and test solutions as described in the regulation. A second stock of test solutions was prepared at a concentration level five times higher than the limits reported in Table 1, which represent the acceptable amount of residual solvents in the final product. System compliance, sensitivity, precision, robustness and linearity were assessed for both water-soluble and water-insoluble pharmaceutical products according to USP <467> method, procedures A, B, C. HS-GC-FID operating parameters, as well as the chromatographic columns, are reported in Table 2. Data was acquired, processed and reported using the Thermo Scientific™ Chromleon™ Chromatography Data System (CDS) software, version 7.2, a platform compliant with Title 21 of the Code of Federal Regulations, Part 11 (Title 21 CFR Part 11). Simplified e-workflows deliver effective data management, ensuring ease of use, data integrity, and traceability. Moreover several ready-made templates are available for the assessment of ICH method validation procedures.

Table 1. Concentration limits in ppm for Class 1, Class 2A and Class 2B residual solvents

Compound Name	Concentration Limit (ppm)	Compound Name	Concentration Limit (ppm)
<b>Class 1</b>		<b>Class 2 A</b>	
1,1-Dichloroethane	8	Methanol	3000
1,1,1-Trichloroethane	1500	Acetonitrile	410
Benzene	2	Dichloromethane	600
Carbon Tetrachloride	4	trans 1,2-Dichloroethane	1870
1,2-Dichloroethane	5	cis 1,2-Dichloroethane	1870
<b>Class 2B</b>		Tetrahydrofuran	720
Hexane	290	Cyclohexane	3880
Nitromethane	50	Methylcyclohexane	1180
Chlorobenzene	60	1,4-Dioxane	380
1,2-Dimethoxyethane	100	Toluene	880
Trichloroethylene	80	Chlorobenzene	360
Pyridine	200	m-Xylene	2170
2-Hexanone	50	p-Xylene	2170
Tetralin	100	o-Xylene	17%

Figure 1. Class 1 system suitability solution peak-to-peak signal-to-noise (S/N) ratios for water-soluble (a) and water-insoluble (b) products.

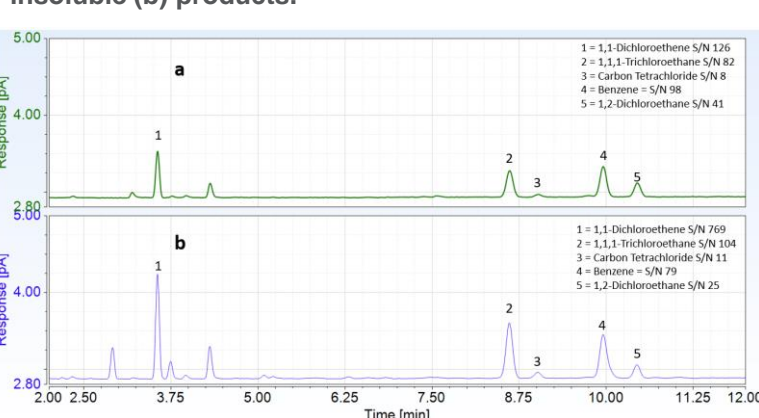
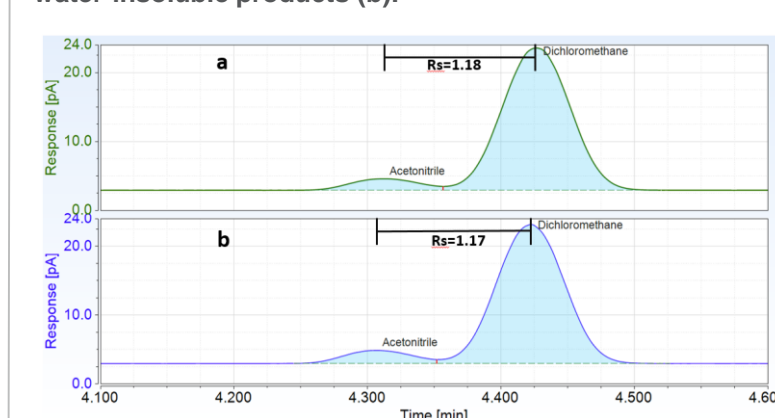


Figure 2. Chromatographic resolution (Rs) between acetonitrile and dichloromethane for water-soluble (a) and water-insoluble products (b).



## RESULTS

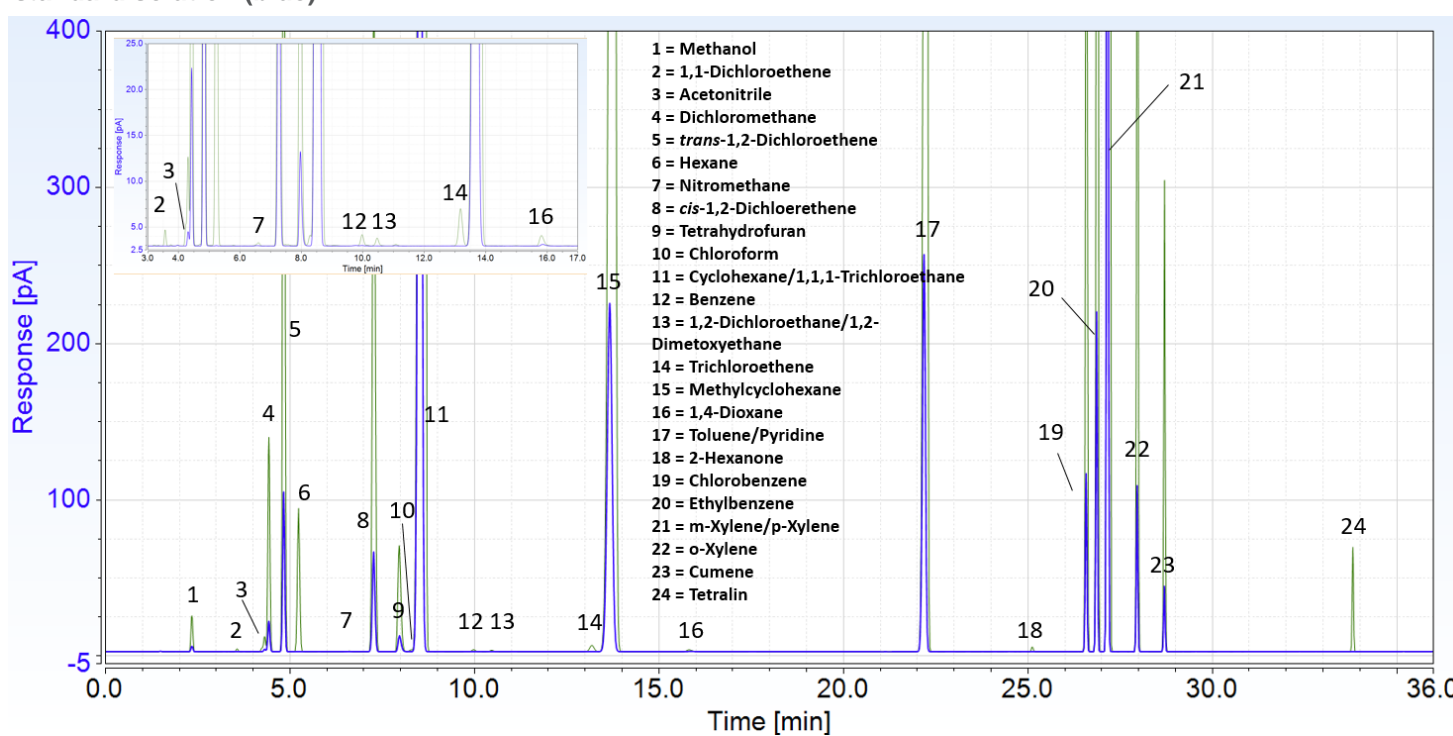
### Procedure A: residual solvent screening and identification

The pharmaceutical products (dispersive aspirin and paracetamol pain relief tablets) and test solutions for water-soluble and water-insoluble pharmaceuticals were prepared as detailed in the USP <467> method and analysed applying the operating conditions reported in Table 2. Peak-to-peak (PP) signal-to-noise ratio (S/N) for 1,1-trichloroethane in Class 1 standard solution was >5:1 and all peaks in Class 1 system suitability showed S/N >3:1 (Figure 1), moreover chromatographic resolution (Rs) between the critical pair, acetonitrile and dichloromethane, was >1 (Figure 2) meeting all the USP <467> method requirements.

### Testing a pharmaceutical product failing procedure A

The pharmaceutical products (dispersive aspirin and paracetamol pain relief tablets) spiked with residual solvents were injected into the chromatographic system. The results were compared to the standard solutions. As peaks found in the spiked samples exceeded the limits reported in Table 1, a compound confirmation step was mandatory as described in the procedure B. As an example, the peak profile obtained for dispersive aspirin spiked solution (green) compared to Class 2A standard solution (blue) is reported in Figure 3.

Figure 3. Comparison between peak profiles obtained for water-soluble spiked test solution (green) and Class 2A standard solution (blue).



Procedure B: peak identity confirmation  
Class 1, Class 1 System Suitability, Class 2A standard solutions, and test solutions for water-soluble and water-insoluble pharmaceuticals were prepared following the USP <467> method and analysed applying the operating conditions reported in Table 2. P/P S/N for benzene in Class 1 standard solution is >5:1 and all peaks in Class 1 system suitability showed S/N >3:1 satisfying the regulation requirements (Figure 4). The critical pair *cis* 1,2-dichloroethane and acetonitrile is baseline resolved with a chromatographic resolution of 3.8 and 3.9 for water-soluble and water-insoluble Class 2A standard solutions, respectively (Figure 5), meeting the required acceptance criteria (Rs ≥ 1.0).

### Testing a pharmaceutical product matching Procedure B confirmation

The peaks identified (procedure A) were confirmed (procedure B) as their responses were higher than the corresponding standards (Figure 6). Therefore, the levels of these residual solvents must be determined (procedure C).

Figure 4. Peak-to-peak signal-to-noise (S/N) ratios for Class 1 system suitability solutions for water-soluble (a) and water-insoluble (b) products.

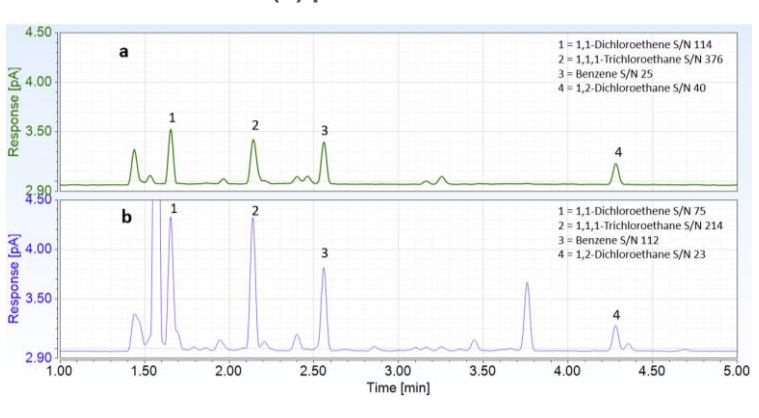
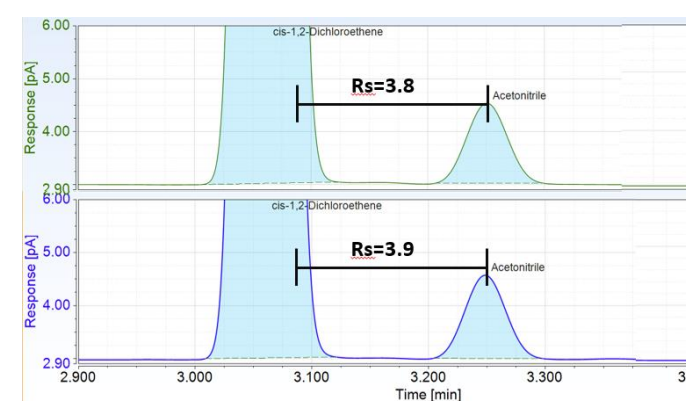


Figure 5. Chromatographic resolution (Rs) between *cis* 1,2-dichloroethane and acetonitrile for water-soluble (a) and water-insoluble products (b).



### Procedure C: quantification

Signal-to-noise (S/N) and chromatographic resolution (Rs) requirements for Class 1, Class 1 System suitability solution, and Class 2A standard solution were the same as described and assessed in procedure A.

### Quantification of the residual solvents in a pharmaceutical product

Class 1, Class 2A, Class 2B standard and test solutions for quantification have been diluted as described by the USP <467> and injected into the chromatographic system. The calculated amount of each residual solvent (in ppm) identified with procedure A and confirmed in procedure B was derived by applying the formula reported in the USP <467> regulation for water-soluble and water-insoluble pharmaceuticals. Calculated concentrations were consistent with the levels used to fortify the samples. As an example, the peak profile for spiked aspirin compared to spiked standard test solution is reported in Figure 7.

Figure 6. Comparison between peak profiles obtained for water-soluble spiked sample solution (green) and Class 2A standard solution (blue).

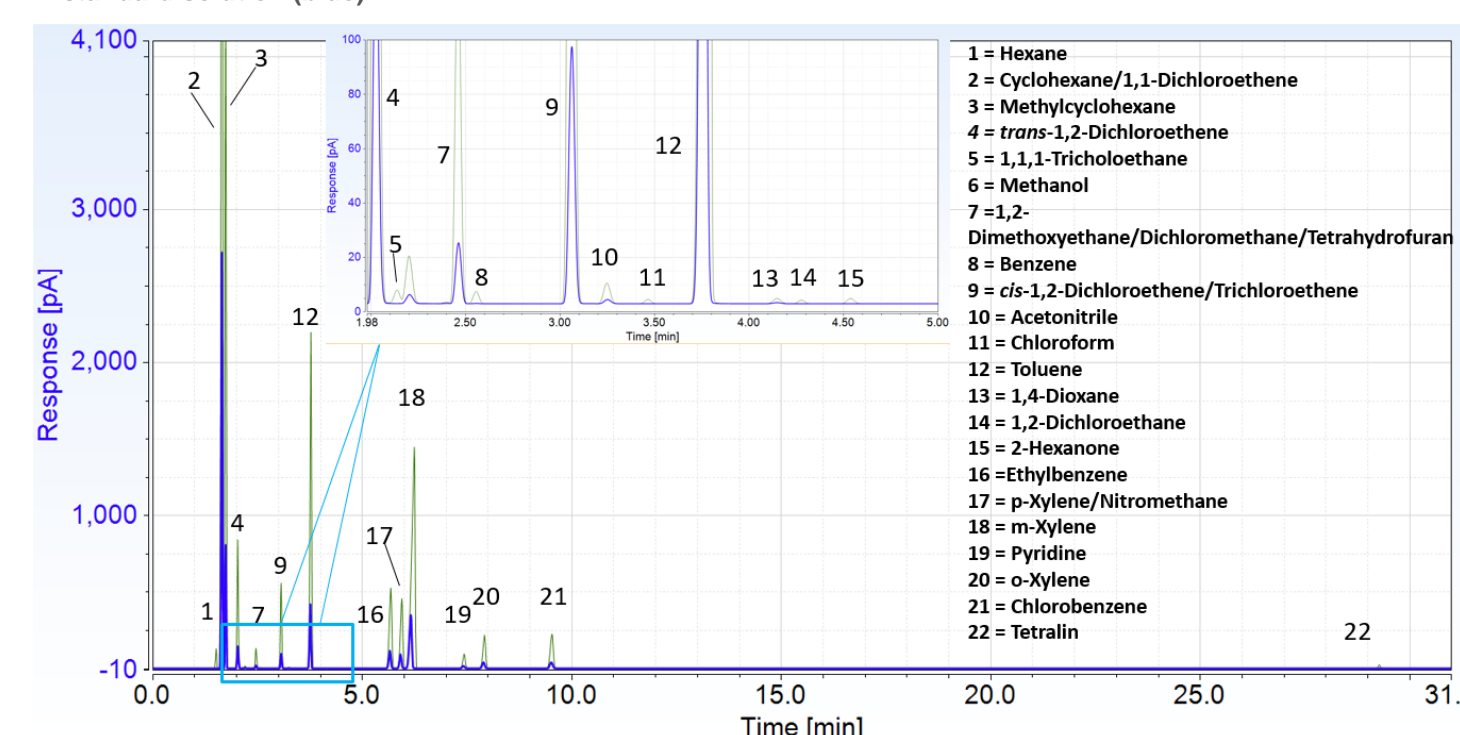
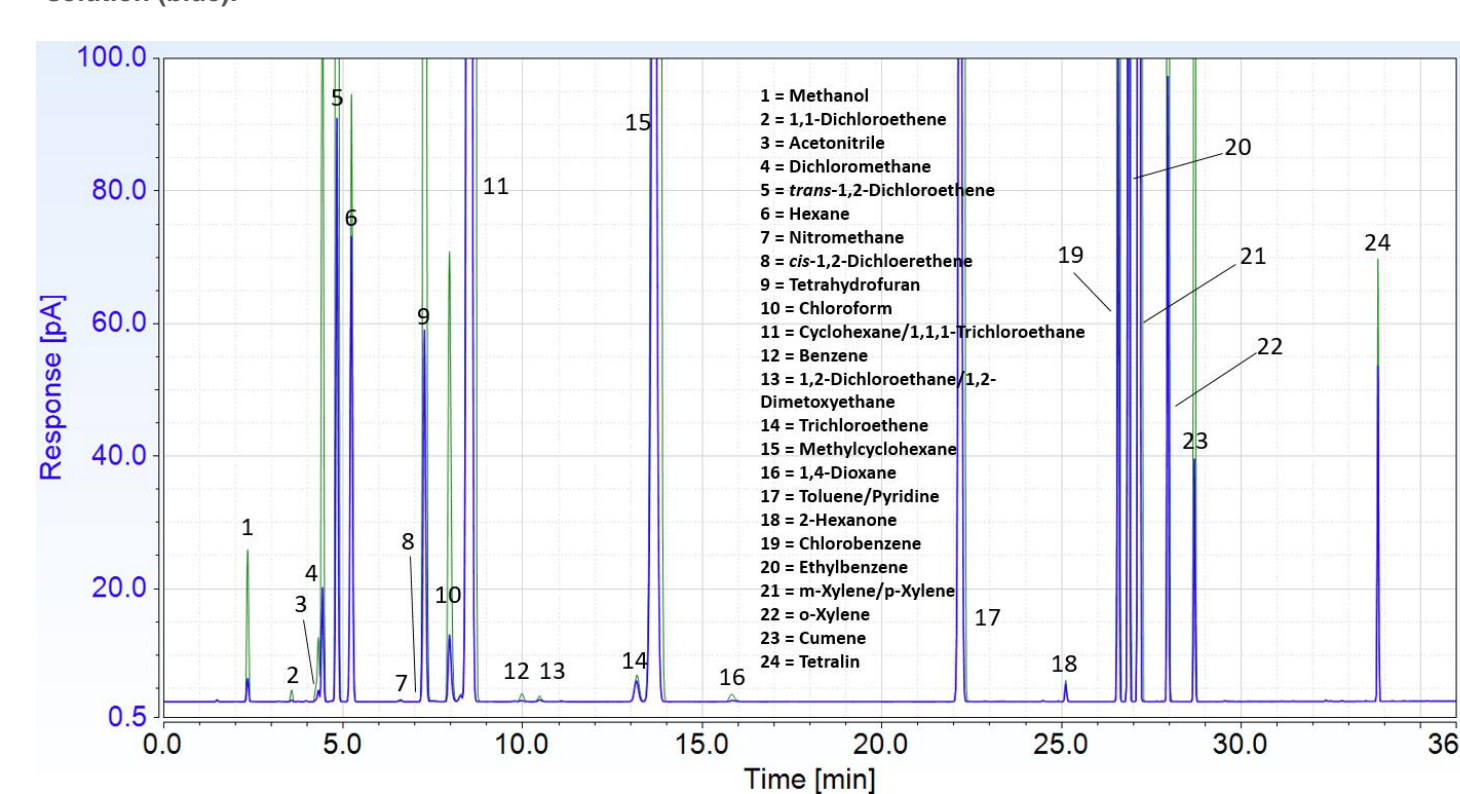


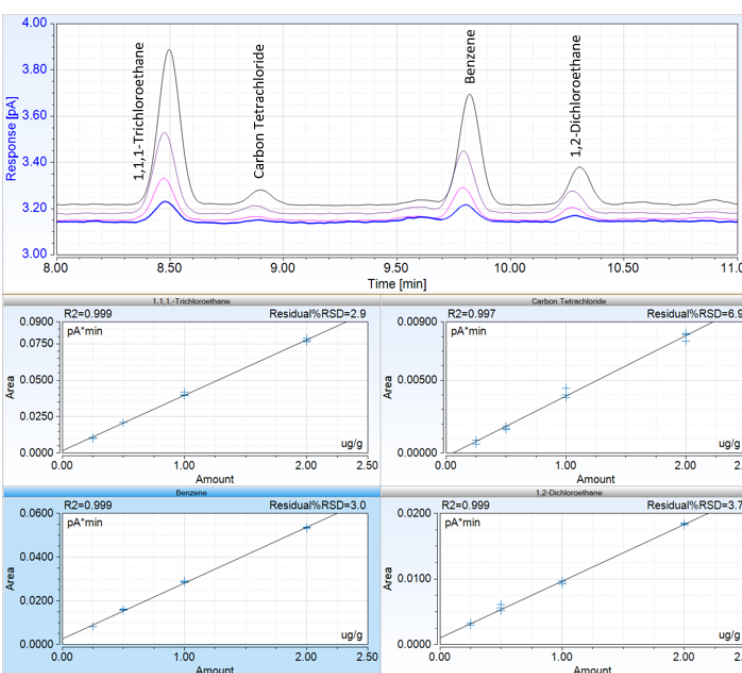
Figure 7. Comparison between peak profiles obtained for spiked aspirin solution (green) and standard test solution (blue).



### Linearity

The standard solutions were diluted to prepare four calibration levels at 12.5%, 25%, 50%, and 100% of the concentration limits reported in the USP <467> method. All residual solvents showed excellent linear responses over the calibration range with an average coefficient of determination: R<sup>2</sup> > 0.999. Moreover, the residual values (measured as % RSD of average response factors) were <4%, confirming a good linear response (Table 3). Each calibration level was prepared and analyzed in triplicate (n=3).

Figure 8. Calibration curves (at 12.5%, 25%, 50% and 100% of the concentration limit) for some of the Class 1 residual solvents.



### Carry-over

Carry-over was assessed analyzing a blank vial after n=9 consecutive injections of pure DMSO (200 µL) and resulted to be <0.0015% (Figure 9). The direct column connection to the valve manifold, the high inert and efficient heated sample path and the effective purging of the loop and the needle contribute to minimize the carry-over effect, especially when high boiling residual solvents are injected.

Figure 9. Carryover <0.0015% after 9 consecutive headspace injections of undiluted DMSO.

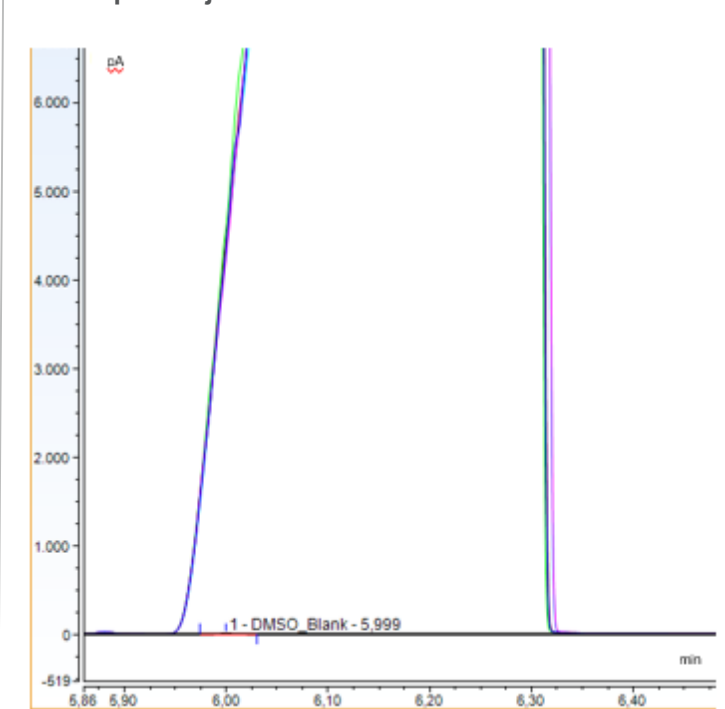


Table 3. Peak area %RSDs obtained from n=18 consecutive injections using water and DMSO as diluents for the concentrated standard solutions, correlation coefficients (R<sup>2</sup>) and relative standard deviation of residuals (%RSD) obtained over four calibration levels at 12.5, 25, 50, and 100% of the concentration limits.

Compound name	%RSD (n=18)		Correlation coefficient (R <sup>2</sup> )	Residuals standard deviation (%RSD)
	Water	DMSO		
1,1-Dichloroethane	1.5	0.7	1.000	2.0
1,1,1-Trichloroethane	1.0	0.8	0.999	2.9
Carbon Tetrachloride	4.9	2.9	0.997	6.9
Benzene	0.8	0.9	0.999	3.0
1,2-Dichloroethane	1.6	1.0	0.999	3.7
Methanol	0.7	1.4	1.000	1.4
Acetonitrile	0.8	1.6	1.000	1.7
Dichloromethane	3.1	0.7	0.998	4.2
trans 1,2-Dichloroethane	4.0	1.2	0.999	2.9
cis 1,2-Dichloroethane	3.4	0.8	0.998	5.0
Tetrahydrofuran	0.9	1.4	1.000	2.2
Cyclohexane	3.6	2.8	0.999	3.0
Methylcyclohexane	3.0	2.4	1.000	2.5
1,4-Dioxane	1.3	1.9	1.000	1.5
Toluene	3.6	0.8	0.997	5.6
Chlorobenzene	3.3	0.7	0.999	2.8
Ethylbenzene	3.4	0.9	0.997	5.3
m-Xylene	3.3	0.9	0.996	6.0
p-Xylene	3.3	0.9	0.996	6.0
o-Xylene	3.1	0.8	0.997	5.6
Hexane	1.2	0.8	0.998	5.8
Nitromethane	2.9	1.5	0.998	4.8
Chloroform	0.9	1.0	0.999	3.0
1,2-Dimethoxyethane	1.4	0.9	0.997	8.4
Trichloroethene	1.9	0.7	0.999	2.9
2-Hexanone	0.6	0.4	1.000	1.3
Tetralin	0.9	0.6	0.999	3.0

Table 5. HS-GC-FID operating parameters

TRACE 1310 GC System Parameters	
Inlet Module and Mode:	SSL split
Split Ratio:	10:1
System Purge Mode, Flow (mL/min):	Constant 5
Carrier Gas, Carrier Mode, Flow (mL/min):	N <sub>2</sub> constant flow, 2.5
Oven Temperature Program:	
Temperature 1 (°C):	40
Hold Time (min):	1
Temperature 2 (°C):	170
Rate (°C/min):	20
FID:	
Temperature (°C):	250
Air Flow (mL/min):	400
H <sub>2</sub> Flow (mL/min):	40
N <sub>2</sub> Flow (mL/min):	40
Acquisition Rate (Hz):	25
Chromatographic Column:	
TG-624 SII MS (P/N 26059-3390):	30 m × 0.32 mm × 1.8 µm
TriPlus 500 HS Autosampler Parameters	
Incubation Temperature (°C):	80
Incubation Time (min):	20
Vial Shaking:	Fast
Vial Pressurization Mode:	Pressure
Vial Pressure (kPa) (Auxiliary Gas Nitrogen):	130
Vial Pressure Equilibration Time (min):	1
Loop Size (mL):	1
Loop/Sample Path Temperature (°C):	80
Loop Filling Pressure (kPa):	73
Loop Equilibration Time (min):	1
Needle Purge Flow Level:	2
Injection Mode:	Standard
Injection Time (min):	1

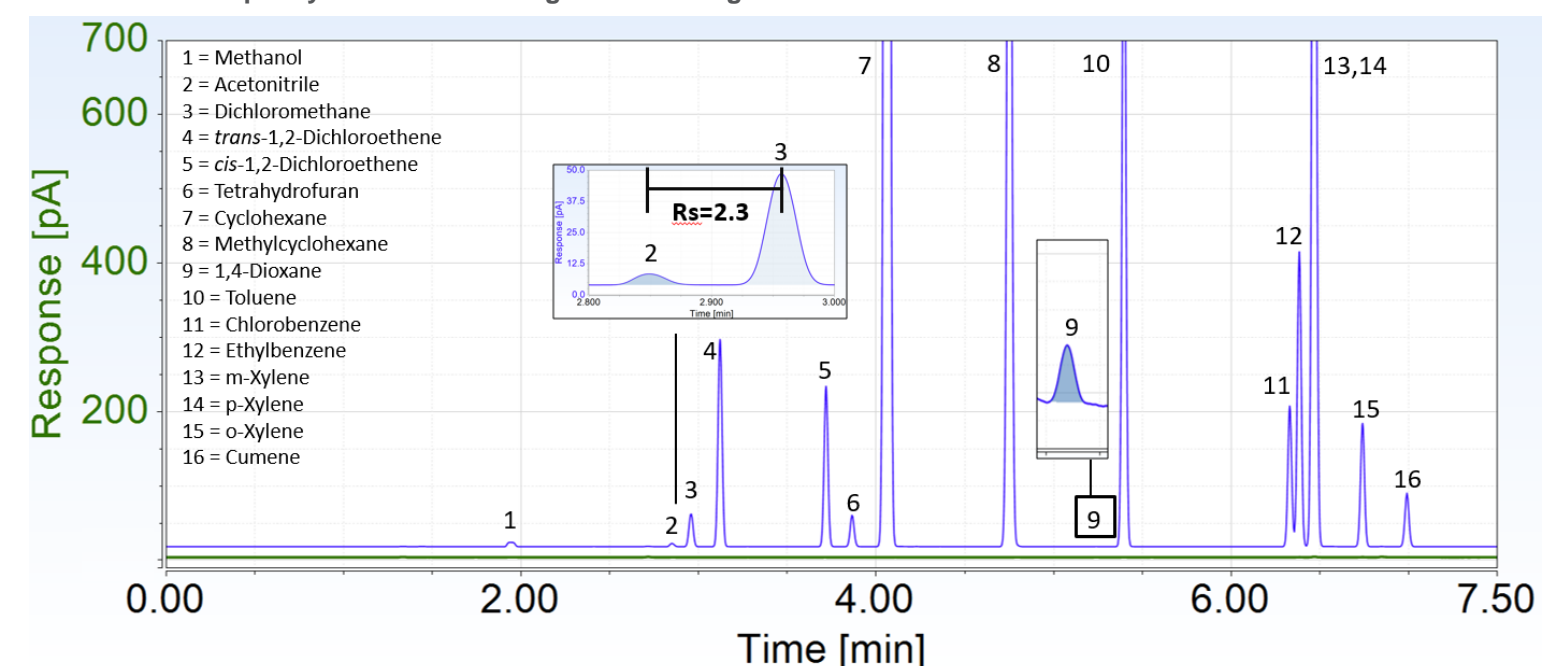
### Repeatability

System repeatability was assessed on n=18 consecutive injections for Class 1, Class 2A, and Class 2B standard solutions. The standard solutions were diluted in water or DMSO according to procedure A for water-soluble and water-insoluble products, respectively. Sample preparation played a critical role for tested apolar solvents with high partition coefficients. As effect of the low affinity for water, %RSDs were higher when concentrated standard solutions were diluted in water with respect to DMSO. Peak area %RSDs obtained for Class 1, Class 2A, and Class 2B residual solvents are reported in Table 3 with average values <3% for all residual solvent classes when water and DMSO were used as diluent.

## SIMPLIFIED AND COST-EFFECTIVE ANALYSIS

The USP General Notices and Requirements<sup>2</sup> allows for the use of alternative methods but they shall be validated as described in the general chapter <1225> Validation of Compendial Procedures. The USP <467> method for Class 2A solvent was modified as an example of how the residual solvent analysis can be improved with a faster and cost-effective alternative method using the TriPlus 500 Headspace autosampler, a USP G43 equivalent capillary column Thermo Scientific™ TraceGOLD™ TG-624 SII MS column and nitrogen as carrier gas. The TG-624 SII MS capillary column allowed for baseline separation of the critical pair acetonitrile/dichloromethane when a fast heating rate was applied, fulfilling the USP <467> resolution requirement (Rs>1) with a short analysis time. Applying a faster GC oven programming rate, the chromatographic separation of Class 2A residual solvents was achieved in less than 8 minutes (Figure 10) as opposed to 60 minutes (with the UPS <467> default conditions), allowing for more than 7 times improvement in analysis speed without compromising chromatographic resolution and method performance. The absolute peak area repeatability was assessed on n=12 consecutive injections of a Class 2A standard solutions at the concentration limits (Table 1). Average peak area %RSD was 1.1% (Table 4), demonstrating a reliable and precise pneumatic control during sampling and injection. Moreover the effective vial shaking allowed for a three times reduction in incubation time without compromising system reliability. HS-GC-FID operating conditions are reported in Table 5.

Figure 10. Chromatographic separation for Class 2A residual solvent can be achieved in less than 8 minutes using a TG-624 SII MS capillary column and nitrogen as carrier gas.



## CONCLUSIONS

- The results presented in this work demonstrate that the new TriPlus 500 HS autosampler in combination with the Trace 1310 GC and FID detector delivers outstanding performance for the analysis of residual solvents in pharmaceutical products meeting or exceeding all USP <467> method requirements.
- The innovative design of the pneumatic control and the flow path inertness ensure excellent repeatability and precision in routine analysis. This was demonstrated by reliable peak area responses obtained (average peak area %RSDs for n=18 consecutive injections was <3%).
- Good linearity (as demonstrated by R<sup>2</sup> and %RSD residual values) was obtained over the calibration range ensuring that the system can be used for routine quantitative assessment of residual solvents in pharmaceutical products.

- The TG-624 SII MS capillary column allowed for fast chromatographic separation of all Class 2A residual solvents in <8 minutes, exceeding the USP <467> chromatographic resolution requirement on the critical pair acetonitrile/dichloromethane (Rs=2.3).
- Sample equilibration was reached in just 20 minutes as a result of the efficient vial spin shaking allowing for a three times reduction in incubation time per sample.

## REFERENCES

- General Chapter USP <467> Organic Volatiles, Chemical Tests, United States Pharmacopoeia, 2012
- General Notices and Requirements USP <38>, United States Pharmacopoeia, 2015
- Thermo Fisher Scientific (2018) Residual Solvent Analysis Application Note 10676
- Thermo Fisher Scientific (2018) Residual Solvent Analysis Technical Note 10679
- Thermo Fisher Scientific (2018) Residual Solvent Analysis Technical Note 10681

## TRADEMARKS/LICENSING

© 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

PO10694-EN 0219S

**ThermoFisher**  
SCIENTIFIC