thermo scientific



Simplified, cost-effective headspace GC method for residual solvents analysis in pharmaceutical products

Authors

Giulia Riccardino and Cristian Cojocariu, Thermo Fisher Scientific, Runcorn, UK

Keywords

USP <467>, residual solvents, pharmaceuticals, valve-and-loop, static headspace, HS, gas chromatography, GC, flame ionization detector, FID, TriPlus 500 Headspace Autosampler

Goal

The aim of this work was to develop a rapid, cost-effective, modified USP <467> HS-GC-FID method for residual solvent determination in pharmaceutical products using the Thermo Scientific[™] TriPlus[™] 500 Headspace Autosampler and nitrogen as carrier gas.

Introduction

Organic solvents are often used in the manufacturing and purification of drug substances but due to their potential toxicity their absence/presence must be verified in the pharmaceutical products to ensure patient safety. The United States Pharmacopeia (USP) method <467>¹ provides detailed procedures for screening, confirmation and quantitation of residual solvents, including sample preparation and analytical conditions.

Gas chromatography (GC) coupled with headspace (HS) sampling technique and flame ionization detection (FID) as detector is the analytical method specified in USP <467> for this application, as most of the target compounds are organic solvents with relatively low boiling points and good thermal stability.

Since a chromatographic resolution (Rs)≥1.0 between the critical pair acetonitrile/dichloromethane is required as a system performance criteria, an initial isothermal condition of the GC column followed by a slow heating rate is recommended by the USP <467> method, in order to avoid peak co-elution and to achieve a baseline peak separation. However, those conditions typically lead to a long analysis time (60 minutes).



The USP General Notices and Requirements² allows for the use of alternative methods but they shall be validated as described in the general chapter <1225> *Validation of Compendial Procedures* and must be shown to give equivalent or better results compared to the USP standard methods.

In this technical note, 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.

Experimental

Sample preparation

USP <467> Class 2A residual solvent solution in dimethylsulfoxide (DMSO) was purchased from Restek[®] (P/N 36012). The stock solution was diluted in DMSO as reported in the USP method <467>, procedure A, water-insoluble articles. HPLC-MS grade water and GC headspace grade DMSO (purity ≥99.9%) were used as diluents. A representative test solution was prepared by diluting locally purchased common pain relief tablets (paracetamol, 500 mg and caffeine, 65 mg) in DMSO, as described in the regulation.

HS-GC-FID analysis

The TriPlus 500 Headspace autosampler was coupled to a Thermo Scientific[™] TRACE[™] 1310 Gas Chromatograph equipped with a Thermo Scientific[™] Instant Connect Split/Splitless (SSL) Injector and a Thermo Scientific[™] Instant Connect Flame Ionization Detector (FID).

Chromatographic separation was achieved on a USP G43 equivalent capillary column Thermo Scientific[™] Trace GOLD[™] TG-624 SiIMS, 30 m × 0.32 mm × 1.8 µm column (P/N 26059-3390) using nitrogen as carrier gas. For this application, nitrogen is a viable alternative to helium (which is expensive and prone to regional shortage). Nitrogen allows for highly efficient GC separations and can easily be produced in the lab at high purity by using a nitrogen generator and making it very cost effective. Full instrumental conditions are given in Table 1.

Data processing

Data was acquired and processed using Thermo Scientific[™] Chromeleon[™] Chromatography Data System (CDS) software, version 7.2, software compliant with the Federal Drug Administration Title 21 Code of Federal Regulations Part 11 (Title 21 CFR Part 11). Table 1. HS-GC-FID analytical parameters for Class 2A residual solvent screening according to the proposed USP <467> method modifications.

TRACE 1310 GC System Parameters		
Inlet Module, Mode	SSL, split	
Split Ratio	10:1	
Septum Purge Mode, Flow	Constant, 5 mL/min	
Carrier Gas, Carrier Mode, Flow	N ₂ , constant flow, 2.5 mL/min	
Oven Temperature Program		
Temperature 1	40 °C	
Hold Time	1 min	
Temperature 2	170 °C	
Rate	20 °C/min	
FID		
Temperature	250 °C	
Air Flow	400 mL/min	
H2 Flow	40 mL/min	
N2 Flow	40 mL/min	
Acquisition Rate	25 Hz	
TriPlus 500 HS Autosampler P	arameters	
Incubation Temperature	80 °C	
Incubation Time	20 min	
Vial Shaking	Fast	
Vial Pressurization Mode	Pressure	
Vial Pressure (Auxiliary Gas Nitrogen)	130 kPa	
Vial Pressure Equilibration Time	1 min	
Loop size	1 mL	
Loop/Sample Path Temperature	80 °C	
Loop Filling Pressure	73 kPa	
Loop Equilibration Time	1 min	
Needle Purge Flow Level	2	
Injection Mode	Standard	
Injection Time	1 min	

Results and discussion

The experiments aimed to improve the analytical method used for residual solvent analysis in pharmaceutical products, and as an example to demonstrate the utility of such method for Class 2A solvents. The ramp is too fast for Class 1. Benzene and 1,2-Dichloroethene co-elute. Furthermore, all the Class 1 RS elute in 10 minutes. The high thermal stability and superior inertness of the TG-624 SiIMS capillary column allowed baseline separation for the critical pair acetonitrile/dichloromethane when a fast heating rate was applied, fulfilling the USP <467> resolution requirement with a short analysis time.

Table 2. Concentration limits (ppm) for Class 2A residual solvents.

Compound	Concentration Limit (ppm)
Class 2A	
Methanol	3000
Acetonitrile	410
Dichloromethane	600
trans 1,2-Dichloroethene	1870
cis 1,2-Dichloroethene	1870
Tetrahydrofuran	720
Cyclohexane	3880
Methycyclohexane	1180
1,4-Dioxane	380
Toluene	890
Chlorobenzene	360
Xylene*	2170
Cumene	70

*Usually 60% m-xylene, 14% p-xylene, 9% o-xylene, 17% ethylbenzene

3

Class 2A standard solution at the concentration limits, as reported in Table 2, was injected into the chromatographic system.

Applying a faster GC oven programming rate, the chromatographic separation of Class 2A residual solvents was achieved in less than 8 minutes (Figure 1) using nitrogen as carrier gas, 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 TG-624 SilMS column allows for excellent chromatographic separation of target compounds, meeting (and significantly exceeding) USP <467> resolution acceptance criteria (Rs≥1.0).

The test solution obtained by diluting common pain relief tablets was injected into the chromatographic system and screened for residual solvent content (Figure 1). No residual solvent peaks could be detected in the pharmaceutical sample, therefore USP <467> requirements for residual solvent content were met (this was in agreement with the same result when using the standard method shown elsewhere.⁴



Figure 1. Chromatographic separation for Class 2A residual solvent can be achieved in less than 8 minutes using a TG-624 SiIMS capillary column and nitrogen as carrier gas. Rs for all residual solvent peaks is ≥1.0, in particular Rs between acetonitrile and dichloromethane met the regulation requirements with a calculated value of 2.3.

Obtaining repeatable data

Vial incubation oven, sample loop and sample path can be set at the same temperature, as the innovative design of the HS-GC direct interface ensures consistent analyte transfer, due to the higher inertness and shorter flow path compared to an external headspace transfer line. Moreover, the precise temperature control of the incubation oven and the effective vial shaking allows for a shorter incubation time (20 minutes) compared to the standard USP <467> method (60 minutes). Importantly, the results precision of target analytes was not affected by speeding up the analysis. This was demonstrated by assessing the absolute peak area repeatability obtained from n=12 consecutive injections of a Class 2A standard solutions at the concentration limits (Table 3). Average peak area %RSD was 1.1%, demonstrating a reliable and precise pneumatic control during sampling and injection.

Table 3. Peak area %RSDs obtained from n=12 consecutive analyses Class 2A residual solvents at the concentration limits described above (Table 2).

Compound	%RSD (n=12)
Methanol	1.3
Acetonitrile	1.2
Dichloromethane	0.8
trans-1,2-Dichloroethene	1.1
cis-1,2-Dichloroethene	0.9
Tetrahydrofuran	1.0
Cyclohexane	1.8
Methycyclohexane	1.5
1,4-Dioxane	1.3
Toluene	1.0
Chlorobenzene	0.8
Ehylbenzene	1.0
<i>m</i> -Xylene/ <i>p</i> -Xylene	1.0
o-Xylene	0.9
Cumene	1.0
Average %RSD	1.1

Conclusions

- Using the experimental conditions described the modified USP <467> method is a real alternative to classical conditions for residual solvent analysis offering the advantages of:
 - a rapid (×7 improvement in analysis speed); cost effective (with nitrogen as carrier gas).
 - high throughput analysis (up to 240 samples analyzed uninterrupted in a single sequence).
- Sample equilibration was reached in just 20 minutes as a result of the efficient vial agitation—a potential three times reduction in incubation time per sample.
- The TG-624 SiIMS 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).
- Reliable pneumatic control during sampling and injection, high inertness of the sample path and consistent analyte transfer through the HS-GC direct interface ensured excellent repeatability of the analyte peak area response, with average peak area %RSD=1.1% from n=12 consecutive vial analyses.
- Chromeleon CDS (compliant with the FDA Title 21 CFR Part 11 requirements) ensures data integrity, traceability, and effective data management allowing for easy and fast data processing, quantitation and reporting.

Taken together, the results from the experiments conducted show that method optimization with nitrogen as carrier gas and TG-624 SilMS capillary column together with Chromeleon CDS represents a viable alternative for routine laboratories.

thermo scientific

References

- 1. General Chapter USP <467> Organic Volatile impurities, Chemical Tests, United States Pharmacopeia, 2012
- 2. General Notices and Requirements USP <38>, United States Pharmacopeia, 2015
- 3. General Chapter USP <621> Chromatography, Physical Tests, United States Pharmacopeia, First Supplement to USP 40-NF 35, 2017
- 4. Thermo Fisher Scientific (2018) Residual Solvent Analysis Application Note: https://assets.thermofisher. com/TFS-Assets/CMD/Application-Notes/an-10676-hsgc-residual-solvents-pharmaceuticals-an10676-en.pdf

Learn More Today

For more information about the TriPlus 500 Headspace Autosampler, to request a demo, or to discuss your GC/GC-MS needs, please contact your Thermo Scientific representative today.

© 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Restek is a registered trademarks of Restek Corp. 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 manners 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 representatives for details. **TN10681-EN 1218M**

