

# All About USP <467> & Residual Solvents

## Regulatory and Application Updates

Manu Grover  
Business Development Manager  
Pharmaceutical Small Molecules Market  
Santa Clara, US

Date: 21<sup>st</sup> Jan. 2025

DE-003843



# Pharmaceutical Impurities

## What

- Unwanted chemicals (either introduced directly using contaminated raw materials or are formed during the mfg. process) that remain with APIs or Drug Product formulations
- Three major categories of pharmaceutical impurities, viz. -
  - Organic & Inorganic Compounds
  - Elemental Impurities
  - **Residual Solvents**

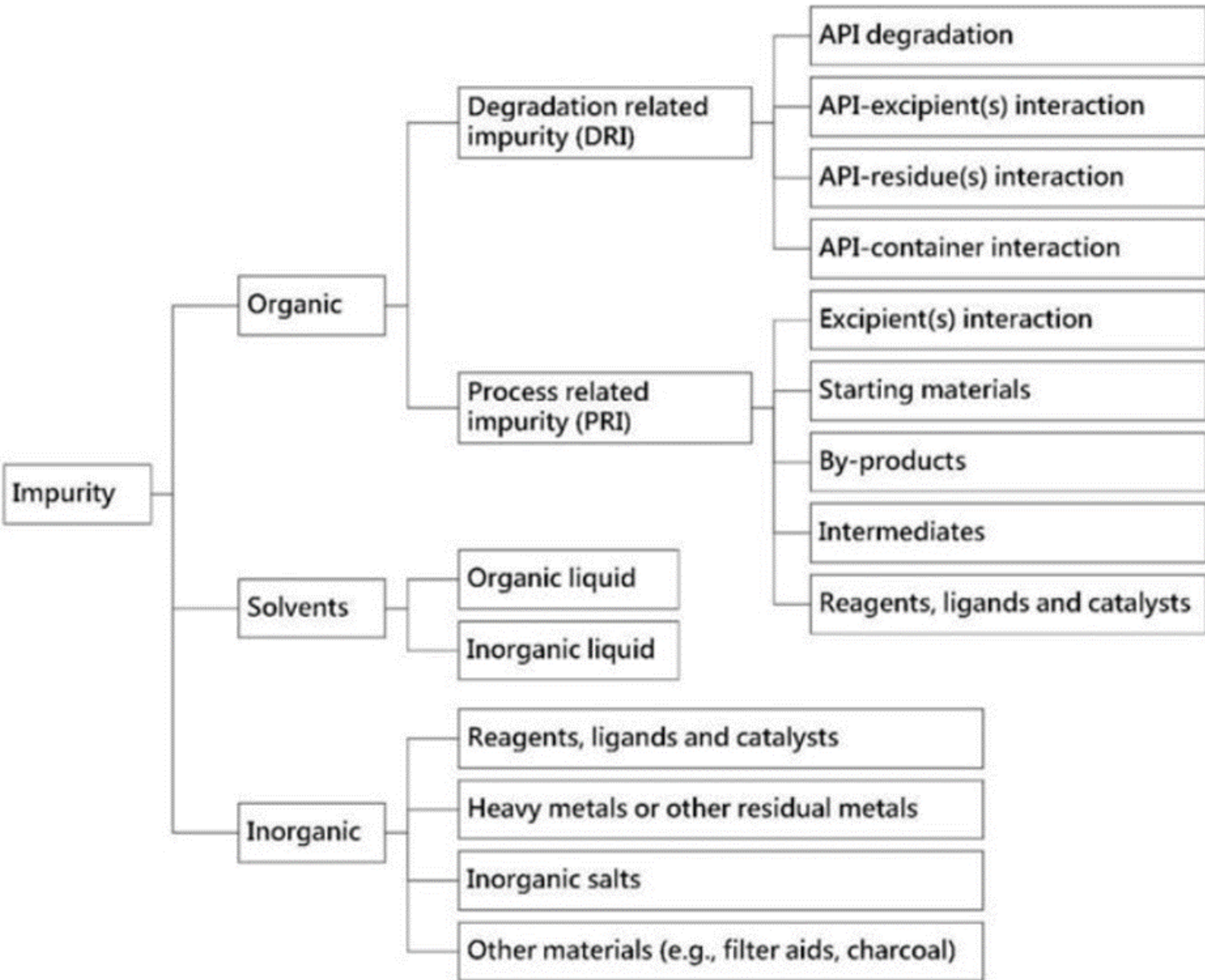
## How

- May arise during synthesis or as a reaction by-product
- Formed as a result of the inherent instability of drug substances
- Added excipients, water
- Interactions with manufacturing equipment and packaging materials and container closure systems (CCSs)

## Why

- The amount of impurity present in drug substance affects the quality and efficacy of the final pharmaceutical product and safety. Therefore, the identification, quantitation, qualification, and control of impurities are a critical part of the drug development process

# Impurity Profiling (Pharmaceuticals)



Ref: (7) (PDF) Determination of Impurities in Pharmaceuticals: Why and How? (researchgate.net)

# What are Residual Solvents?

**Residual solvents** – Organic volatile chemicals used or produced in manufacturing of API's, excipients or formulations.

Residual solvents can arise from various sources:

- Synthesis or Reaction byproduct
- Inherent instability exhibited by some drug substances
- Excipients and water used in the manufacturing process
- Interactions with manufacturing equipment and packaging materials, including container closure systems.

The primary guiding regulations for residual solvents include [USP <467>](#), which aligns closely with [ICH Q3C\(R8\) guidelines](#).

*'Since there is **no therapeutic benefit** from residual solvents, these should be removed to extent possible to meet Pdt Specs, GMP or other requirements.'*

The document classifies residual solvents into 3 categories:

- **Class 1:** Solvents to be avoided, *including known human carcinogens, strongly suspected human carcinogens, and solvents posing environmental hazards.*
- **Class 2:** Solvents to be limited, *comprising nongenotoxic animal carcinogens or agents causing other irreversible toxicities such as neurotoxicity or teratogenicity. This class also includes solvents suspected of causing significant but reversible toxicities.*
- **Class 3:** Solvents with low toxic potential, *having no health-based exposure limit.*

# Residual Solvent Analysis

USP <467> and ICH Q3C (R5 and R8) compliance for residual solvent analysis

Organic volatile chemicals used in the preparation of APIs, excipients & drug products

ICH

USP

Require Chromatographic Analysis



- Solvents to be avoided
- Know or strongly perceived human carcinogens

Class 1



- Solvents to be limited
- Non-genotoxic carcinogens or possible neurotoxic or teratogenic agents

Class 2



- Solvents with low toxic potential
- No health-based exposure limit is needed

Class 3





# Residual Solvents limits & Control Strategies

**Permitted Limits** – Method used to establish Permitted Daily Exposure (PDE) for residual solvents is derived from no-observed effect level (NOEL) or lowest-observed effect level (LOEL) in most relevant animal study (Toxicity Data).

*‘Class 1 solvents should be avoided, Meanwhile limits for Class 2 solvents are based on the toxicological permitted daily exposure (PDE), whereas Class 3 solvents are considered less toxic, and control to 50 mg/day or less for each of these solvents is acceptable without justification.’*

Options to describe limits of solvents:

**Option-1:** Concentration (ppm) =  $1000 \times \text{PDE} / \text{Dose}$

**Option-2:** Adding the amount of residual solvents present in each component of drug product. The SUM of amounts of solvent per Day should be less than PDE.

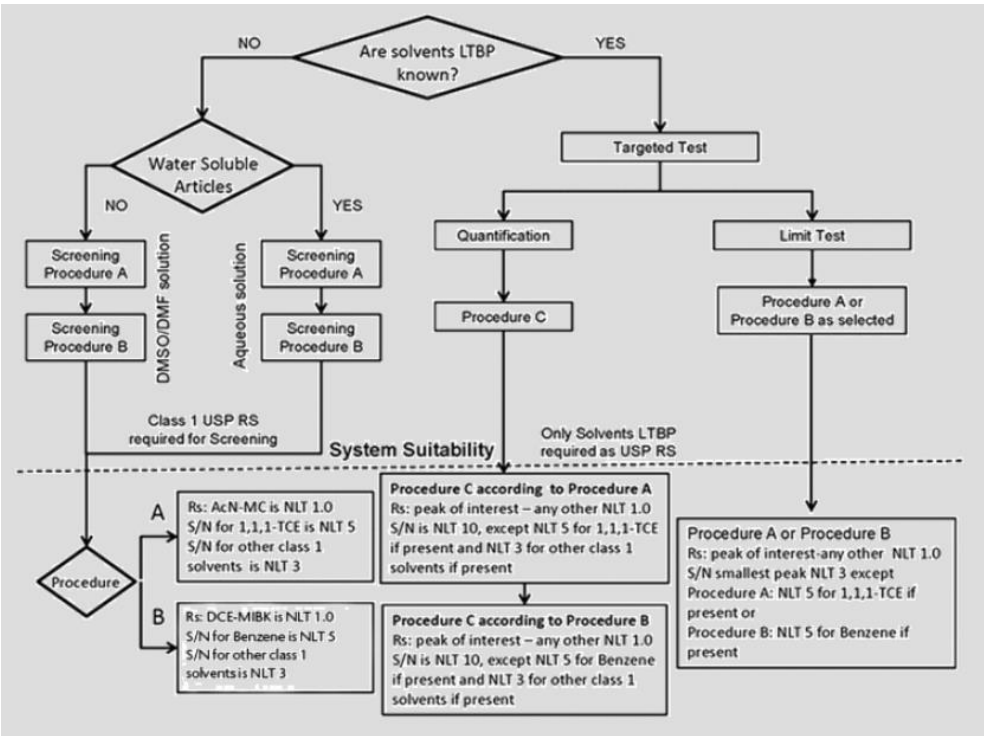
## ***Guidelines for Residual Solvents (CPMP/ICH/283/95)***

- When a Class 1 solvent might be present in another solvent (e.g. Acetone or Toluene containing Benzene), routine testing is not required when – The limit applied to original solvent is such that class 1 solvent will be present in API at levels below guidelines limits.
- It is proved (with Validated method) that class 1 solvent is not more than 30% of spec limits, in intermediates or final pdt.
- Supporting data – 6 consecutive pilot scale or 3 consecutive Industry scale batches.
- When a Class 2 is LTBP, they should be routinely controlled in intermediate or final pdt. e.g.
  - LTBP in last step of synthesis it should be routinely controlled in Final API/ Pdt.
  - LTBP prior to last step, then need not to be reported in Drug Specs if levels are NMT 10% of acceptable limits (e.g. Acetonitrile 41 ppm)

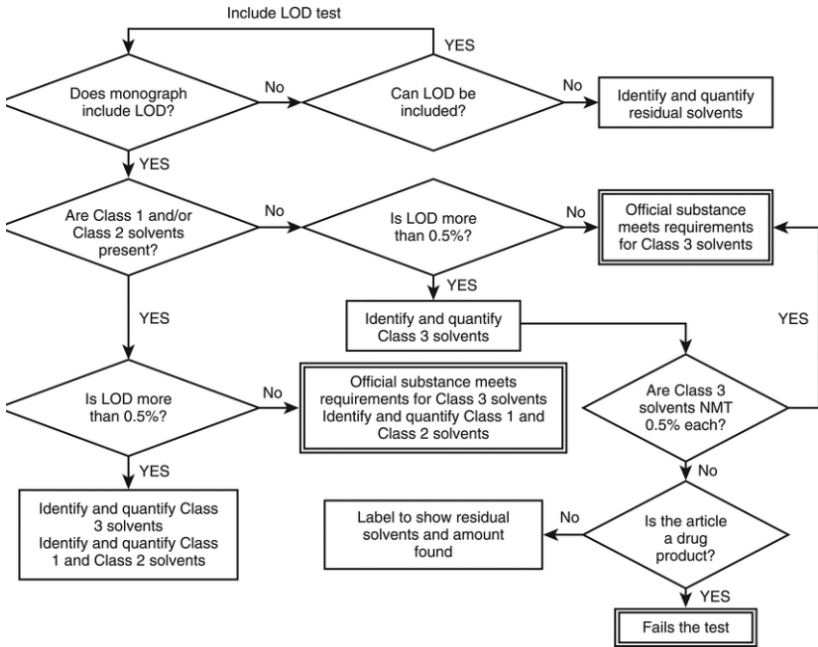
# Residual Solvents limits & Control Strategies

<467> describes Analytical procedures (Procedures A, B, and C) for evaluating the levels of all Class 1 and the majority of Class 2 residual solvents. For each test matrix, verification is needed to demonstrate the reliability of the compendial procedure, as described in <1467>.

- When the solvents **LTBP (likely to be present) are not known**, use the screening tests of Procedure A and B, as required.
- When the information about solvents **LTBP in the material is known**, the system suitability requirements only need to be demonstrated for the solvents expected to be present.



USP<467> Figure 4. System suitability requirements based on the method of choice and previous knowledge about solvents LTBP.



USP<467> Figure 5. Decision tree for the control of Class 3 residual solvents

# Limits of Residual Solvents

TABLE 1: CLASS 1 RESIDUAL SOLVENTS<sup>1</sup>

| RESIDUAL SOLVENT      | CONCENTRATION LIMIT (PPM) | TOXICOLOGICAL CONCERN          |
|-----------------------|---------------------------|--------------------------------|
| Benzene               | 2                         | Carcinogen                     |
| Carbon Tetrachloride  | 4                         | Toxic and Environmental Hazard |
| 1,2-Dichloroethane    | 5                         | Toxic                          |
| 1,1-Dichloroethene    | 8                         | Toxic                          |
| 1,1,1-Trichloroethane | 1500                      | Environmental Hazard           |

TABLE 3: CLASS 3 RESIDUAL SOLVENTS -

| RESIDUAL SOLVENT       |
|------------------------|
| Acetic acid            |
| Acetone                |
| Anisole                |
| 1-Butanol              |
| 2-Butanol              |
| Butyl acetate          |
| Tert-Butylmethyl ether |
| Cumene                 |
| Dimethyl sulfoxide     |
| Ethanol                |
| Ethyl acetate          |
| Ethyl ether            |
| Ethyl formate          |
| Formic acid            |
| Formic acid            |
| Ethyl formate          |
| Ethyl ether            |
| Ethyl acetate          |

|                      |
|----------------------|
| Heptane              |
| Isobutyl acetate     |
| Isopropyl acetate    |
| Methyl acetate       |
| 3-Methyl-1-Butanol   |
| Methylethylketone    |
| Methylisobutylketone |
| 2-Methyl-1-propanol  |
| Pentane              |
| 1-Pentanol           |
| 1-Propanol           |
| 2-propanol           |
| Propyl acetate       |

TABLE 2: CLASS 2 RESIDUAL SOLVENTS, WHERE PDE IS THE PERMITTED DAILY EXPOSURE<sup>2</sup>

| RESIDUAL SOLVENT      | PDE (MG/DAY) | CONCENTRATION LIMIT (PPM) |
|-----------------------|--------------|---------------------------|
| Acetonitrile          | 4.1          | 410                       |
| Chlorobenzene         | 3.6          | 360                       |
| Chloroform            | 0.6          | 60                        |
| Cyclohexane           | 38.8         | 3880                      |
| 1,2-Dichloroethene    | 18.7         | 1870                      |
| 1,2-Dimethoxyethane   | 1.0          | 100                       |
| N,N-Dimethylacetamide | 10.9         | 1090                      |
| N,N-Dimethylformamide | 8.8          | 880                       |
| 1,4-Dioxane           | 3.8          | 380                       |
| 2-Ethoxyethanol       | 1.6          | 160                       |
| Ethylene glycol       | 6.2          | 620                       |
| Formamide             | 2.2          | 220                       |
| Hexane                | 2.9          | 290                       |
| Hexane                | 5.8          | 580                       |
| Formamide             | 5.5          | 550                       |
| Ethylene glycol       | 8.5          | 850                       |
| 2-Ethoxyethanol       | 1.6          | 160                       |

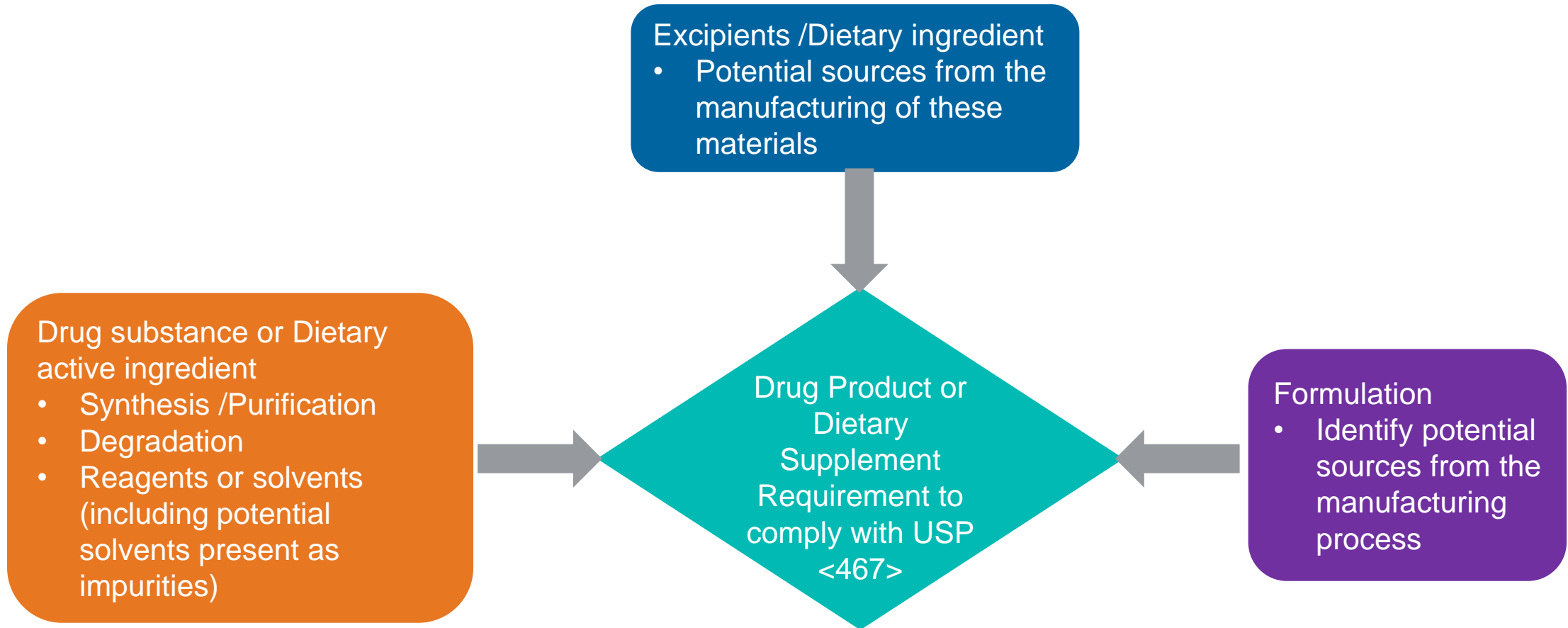
|                     |      |      |
|---------------------|------|------|
| Methanol            | 30.0 | 3000 |
| 2-Methoxyethanol    | 0.5  | 50   |
| Methylbutylketone   | 0.5  | 50   |
| Methylcyclohexane   | 11.8 | 1180 |
| Methylene chloride  | 6.0  | 600  |
| N-Methylpyrrolidone | 5.3  | 530  |
| Nitromethane        | 0.5  | 50   |
| Pyridine            | 2.0  | 200  |
| Sulfolane           | 1.6  | 160  |
| Tetrahydrofuran     | 7.2  | 720  |
| Tetralin            | 1.0  | 100  |
| Toluene             | 8.9  | 890  |
| Trichloroethylene   | 0.8  | 80   |
| Xylene              | 21.7 | 2170 |

Residual solvents should be limited in drug substances, excipients, dietary ingredients, and official products because of the inherent toxicities of these residual solvents. - [467 RESIDUAL SOLVENTS \(uspnf.com\)](https://www.uspnf.com/usp/467/usp467-residual-solvents)



# Potential Sources of Residual Solvents to be Considered

In pharmaceutical drug products and dietary supplements



*[For more details: refer to 467 RESIDUAL SOLVENTS \(uspnf.com\)](https://www.uspnf.com)*

# Changes to Residual Solvents

## USP <467> – Official December 2020

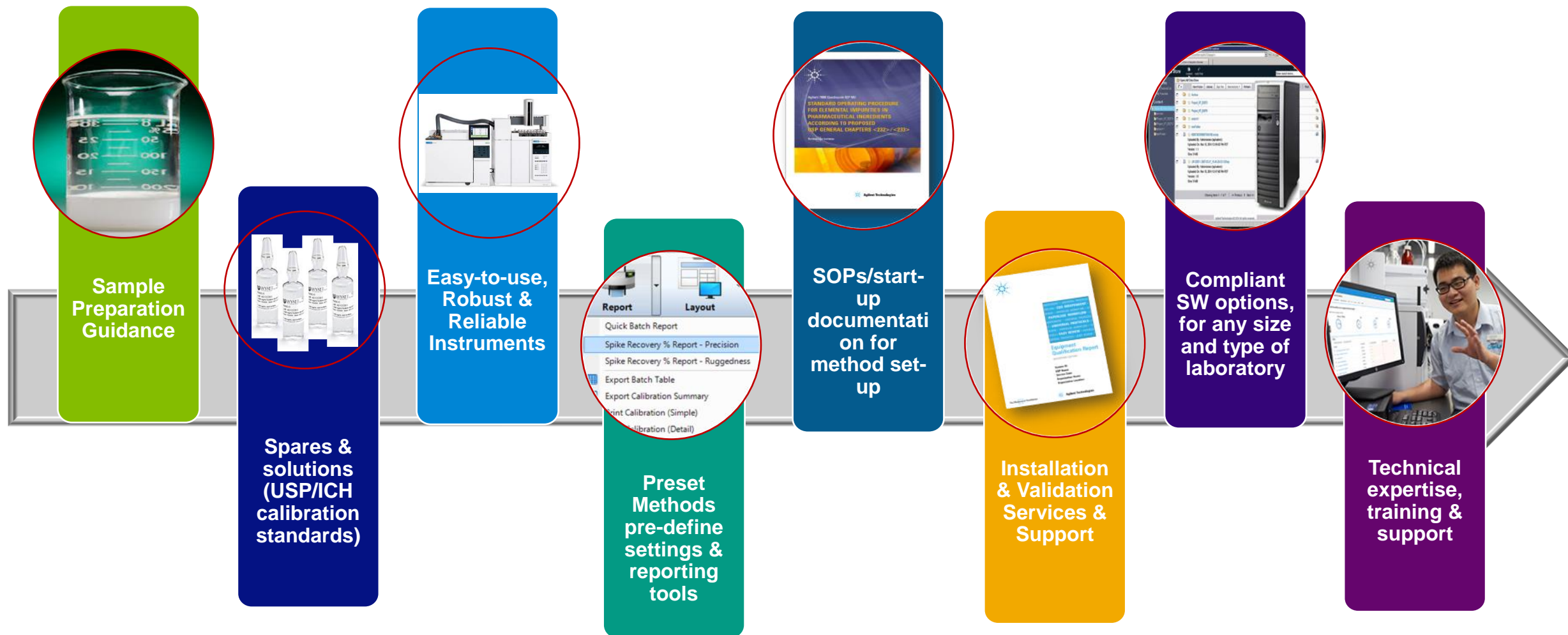
- Methylisobutylketone added to Class 2 list
- Methylisobutylketone used for system suitability resolution

## Three new solvents to be added in next draft

- 2-Methyltetrahydrofuran – Class 3
- Cyclopentyl methyl ether – Class 2
- Tert-butyl alcohol – Class 2



# Agilent Innovative Solutions for Residual Solvent Analysis



# A Flexible Portfolio of Products

Something for everyone!



Ultimate productivity & usability, especially for MS

Intuvo

Quality  
Smart  
Connected

8860

8890



Core routine  
applicability



Advanced  
flexibility &  
expandability

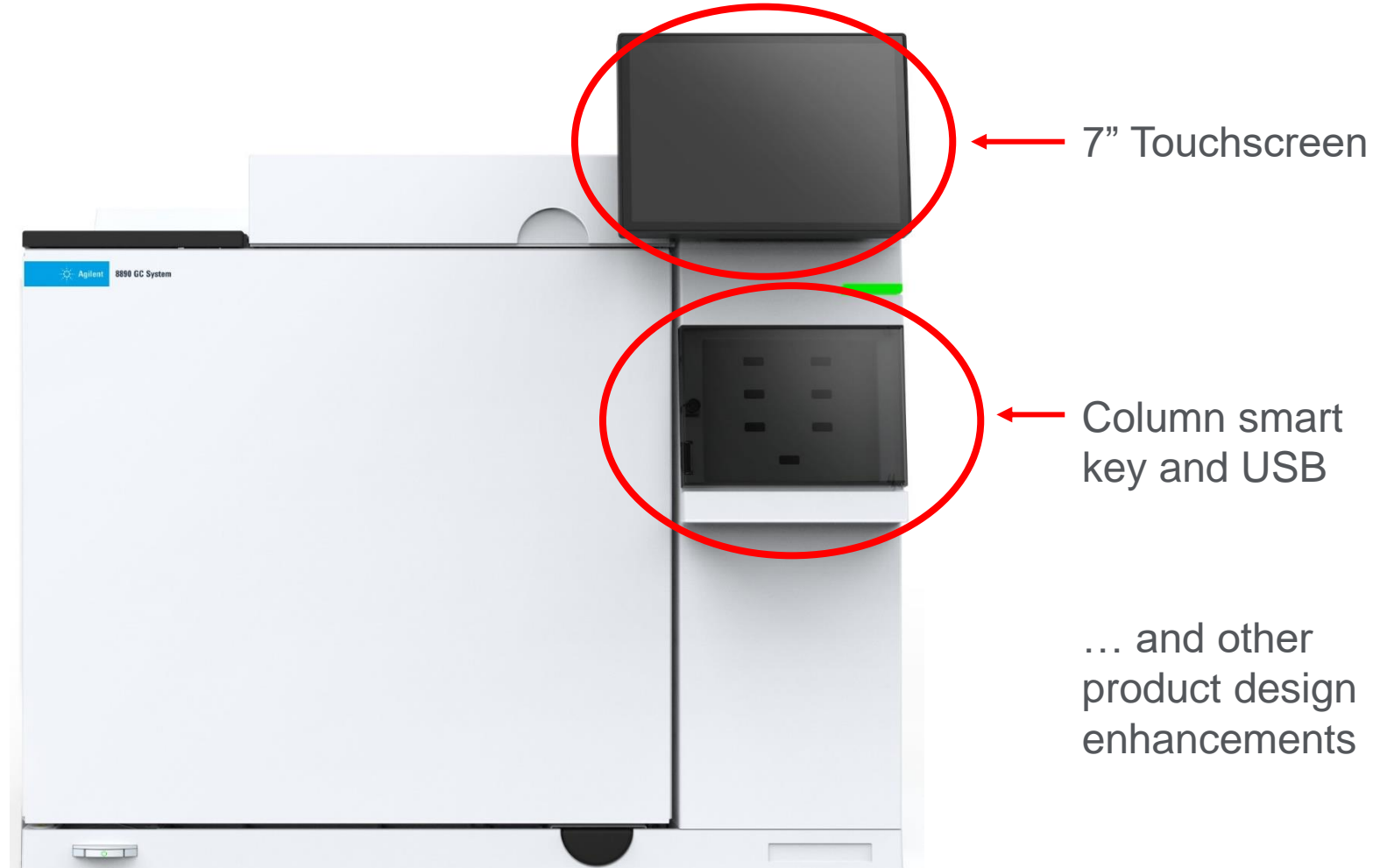
# 8890 Enhancements

## GUI/Display/Touchscreen

- “At-GC”
  - Enhanced Diagnostics
  - Chromatographic Attributes
  - Maintenance
  - Instrument Status
  - Extensive User Info
- Capacitive touch screen

## On-Board CPU → Browser UI

- GC info available via PC, Tablet, Mobile





# Agilent's GC & GC/MSD for Residual Solvent Analysis

Routine Analysis

Rapid High Throughput

Sensitivity

Intuvo 9000/8697HSS



8890GC/7697HSS



8890GC/5977MSD



All systems are equally capable of delivering all USP/ICH compliance requirements

# Agilent's Residual Solvents Workflow

Which GC is right for my analysis?



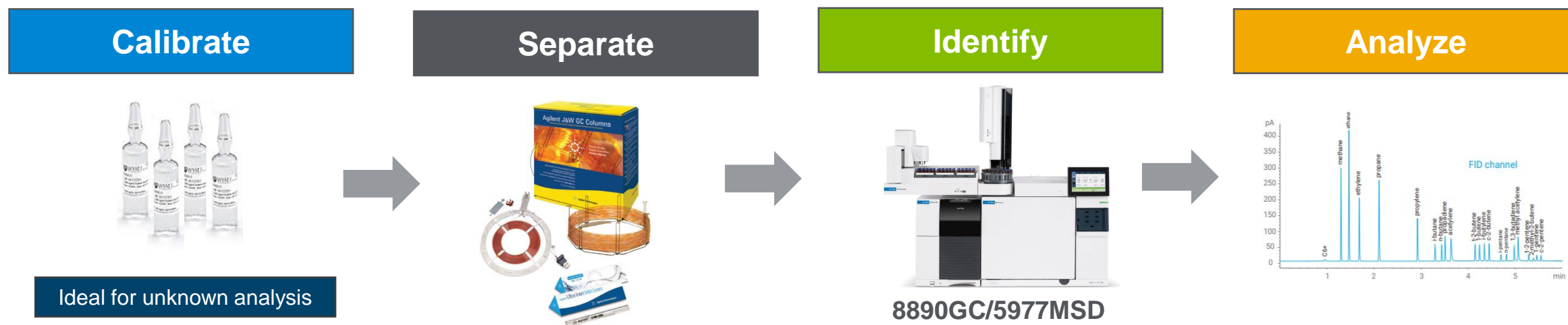
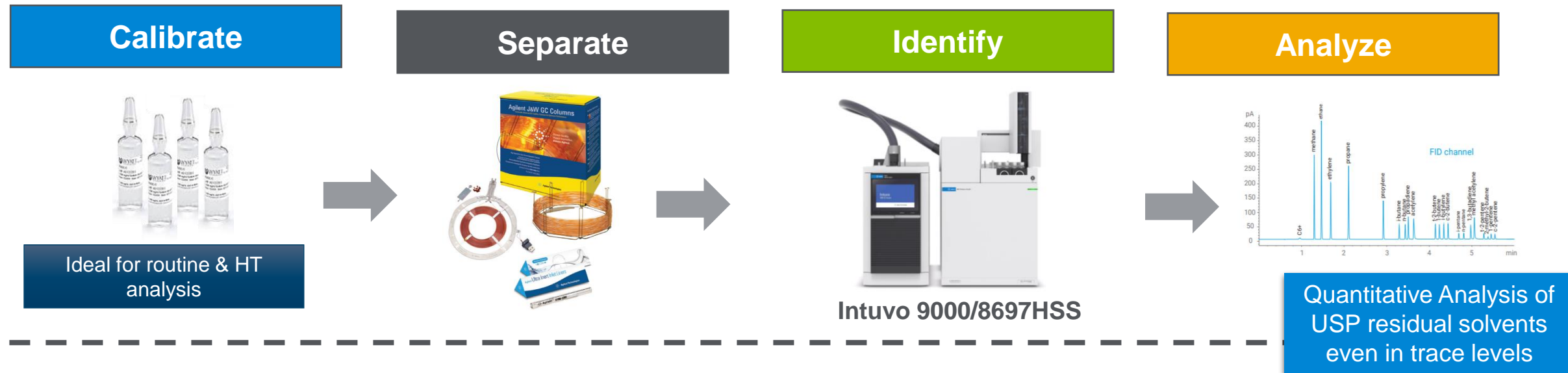
Enhanced detection and identification analysis options of analytes in a single mix -

- Splitting to two columns for dual detector analysis
- Agilent GC coupled to GC/MS System

| GC System / Attributes | Intuvo 9000 | 8890 |
|------------------------|-------------|------|
| Performance            | ✓✓          | ✓    |
| Detector (FID and MS)  | ✓✓          | ✓✓   |
| Footprint              | ✓✓          | ✓    |
| Compliant SW           | ✓✓          | ✓✓   |
| Non-Agilent CDS        | ✓✓          | ✓✓   |
| Direct Replacement     | ✓           | ✓✓   |
| Revalidate SOP         | ✓           | ✓✓   |

# Agilent GC/FID and GC/MS Workflow for Residual Solvent Analysis

Confidently identify and quantify residual solvent in APIs and drug products



# USP <467> Residual Solvent Requirements

USP <467> specifies a single-column analysis

- A secondary analysis is performed if the solvent is found above limit detection.
- An Intuvo 9000 GC configured with an inlet split to two columns and two FIDs can perform both analyses in a single run.



USP 40, general chapter USP <467> Residual Solvents <https://hmc.usp.org/sites/default/files/documents/HMC/GCs-Pdfs/c467.pdf>

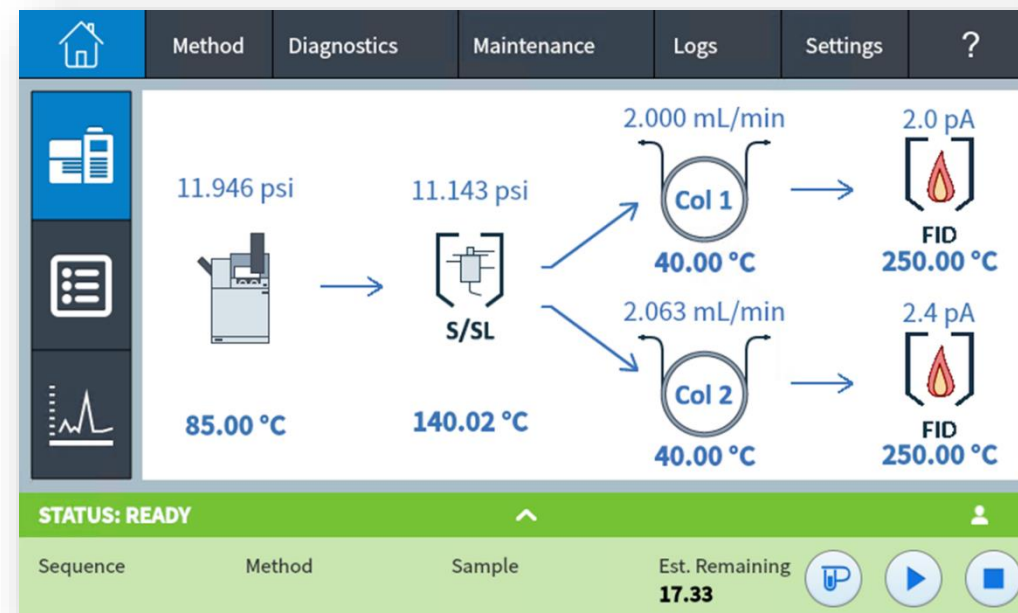
# USP <467> Columns and Performance Requirements

## Procedure A – Initial identification and limit test

- DB-Select 624 UI
- s/n of 1,1,1-trichloroethane > 5
- s/n of all Class 1 solvents > 3
- Resolution of acetonitrile and methylene chloride > 1

## Procedure B – Secondary analysis for confirmation

- DB-Wax Ultra Inert
- s/n of benzene > 5
- s/n of all Class 1 solvents > 3
- Resolution of **methylisobutylketone** and cis-dichloroethene > 1

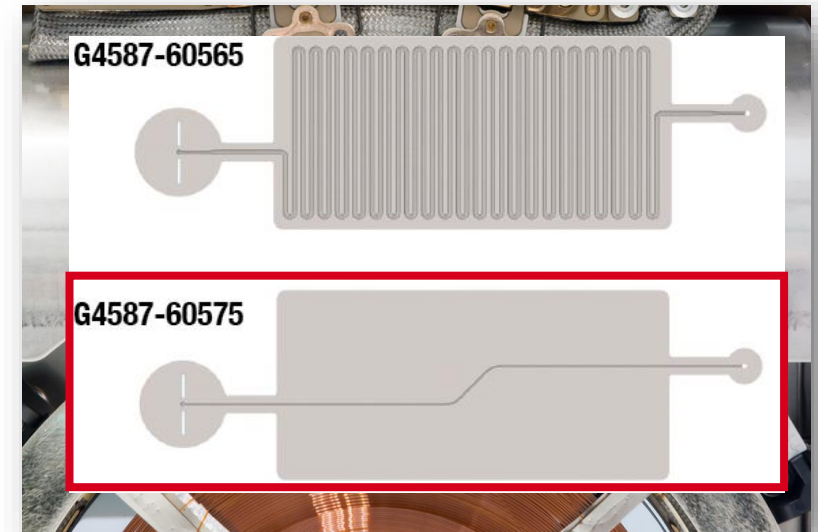
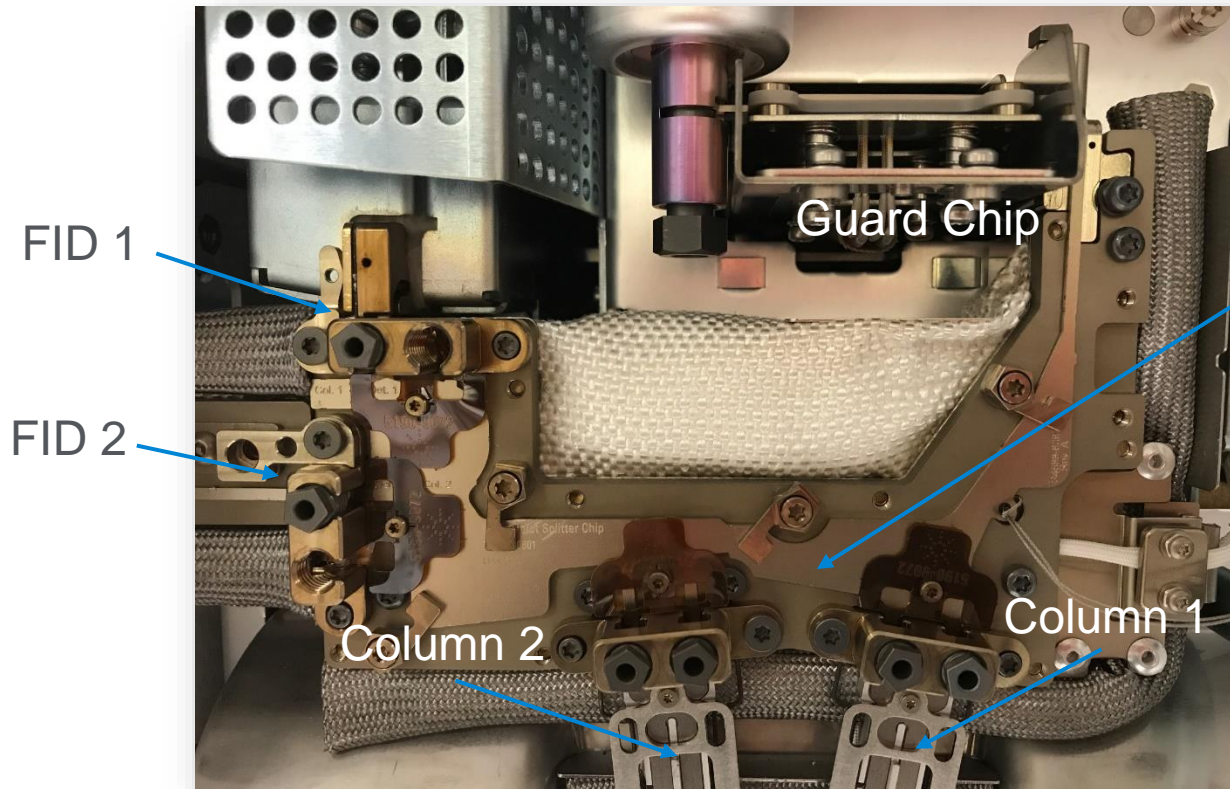




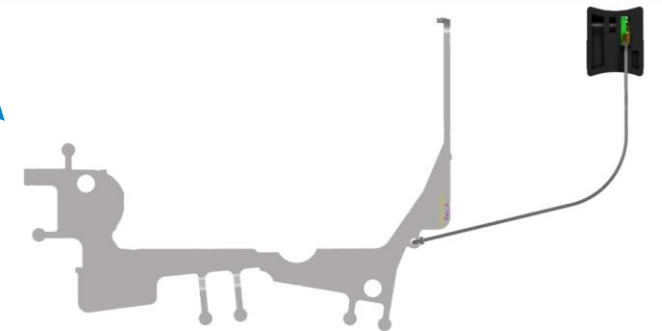
# Configuration

- Headspace to SSL
- Inlet splitter flow chip – equally splits to two columns
- Two columns to two FID detectors

Inlet

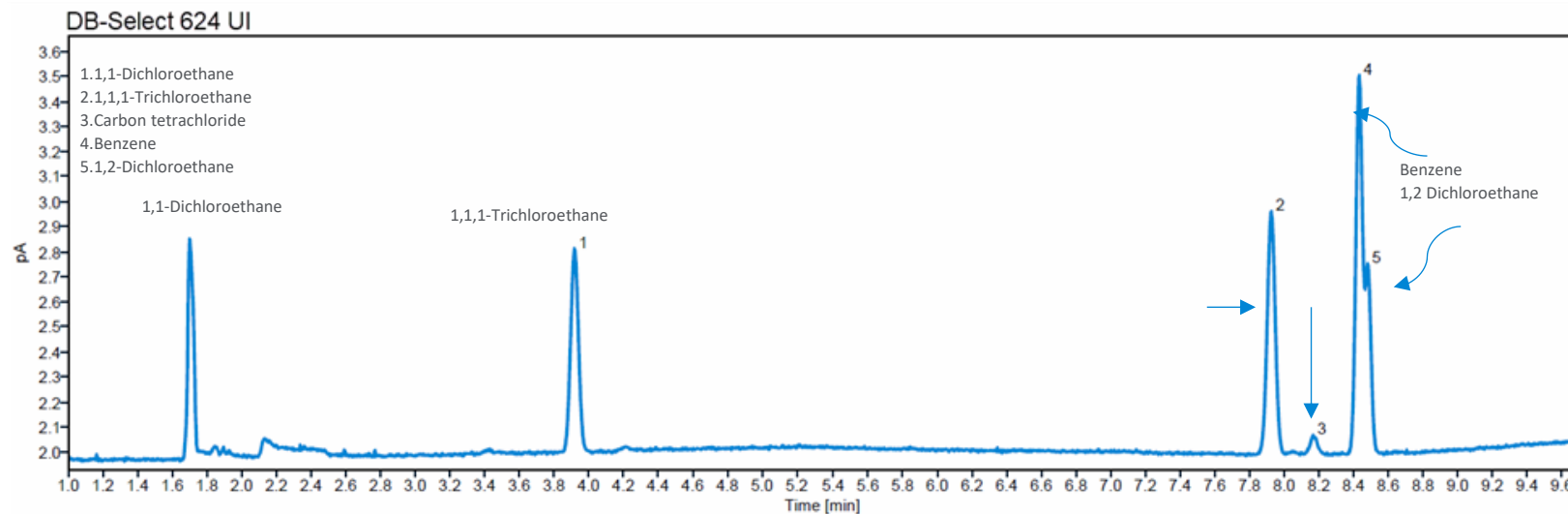


Inlet splitter flow chip

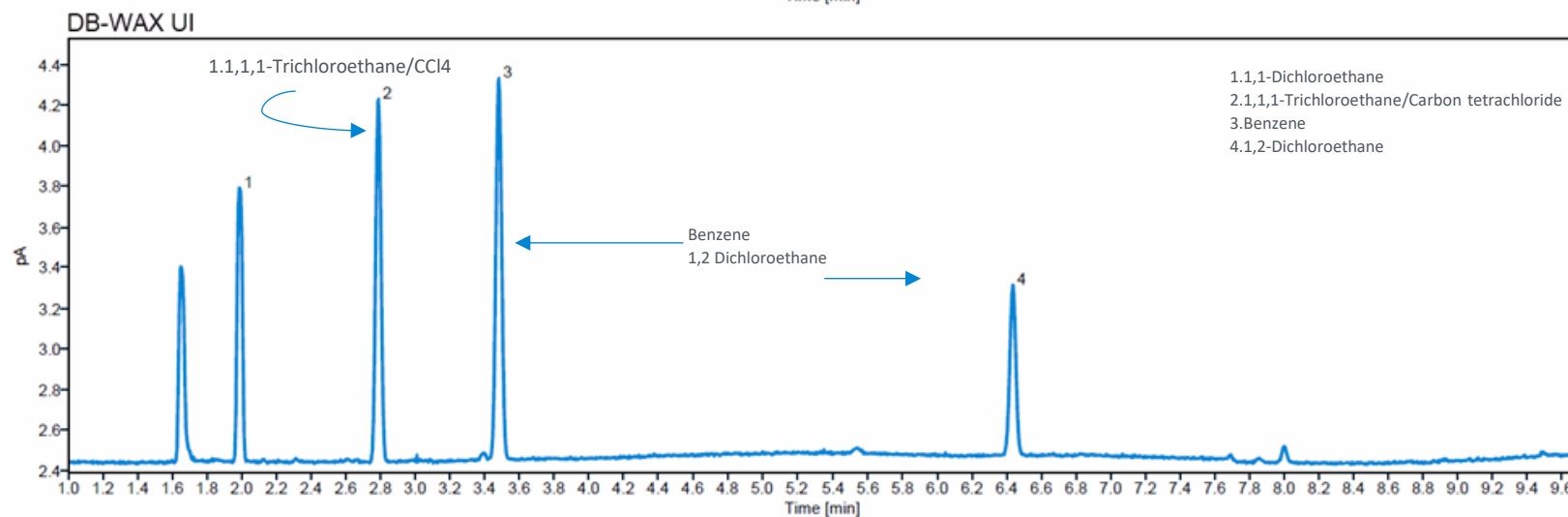


# Class 1, DB-Select 624 UI & DB-WAX Ultra Inert

Signal to Noise  
requirements  
are met

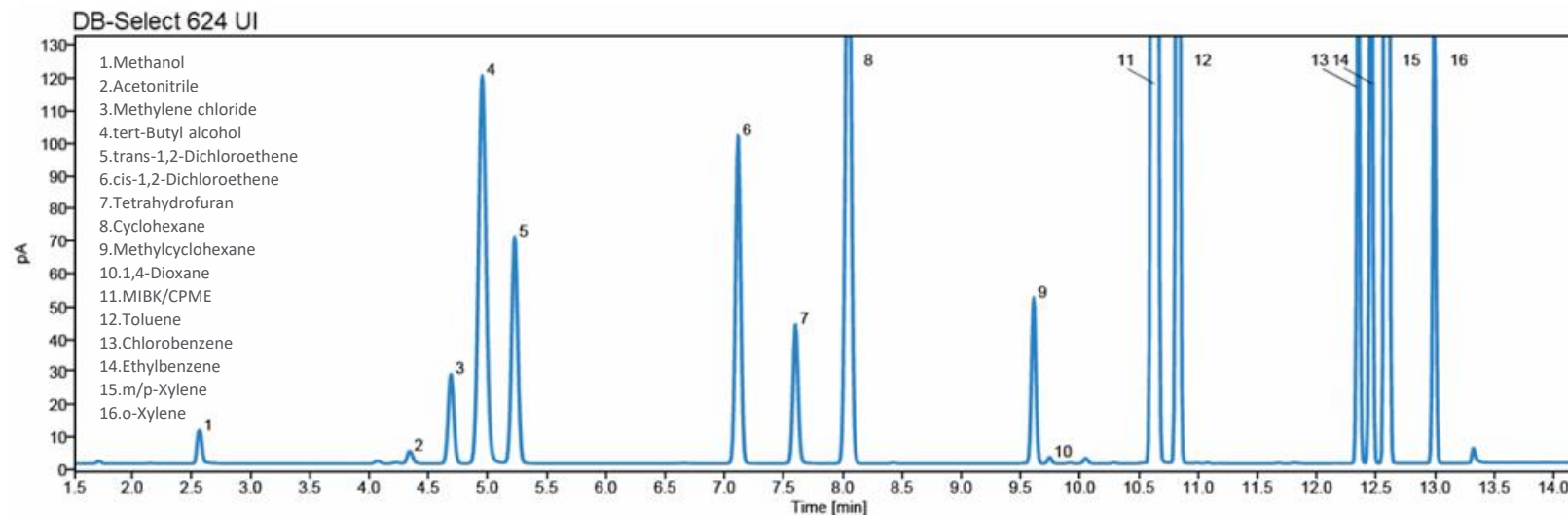


Signal to Noise  
requirements  
are met

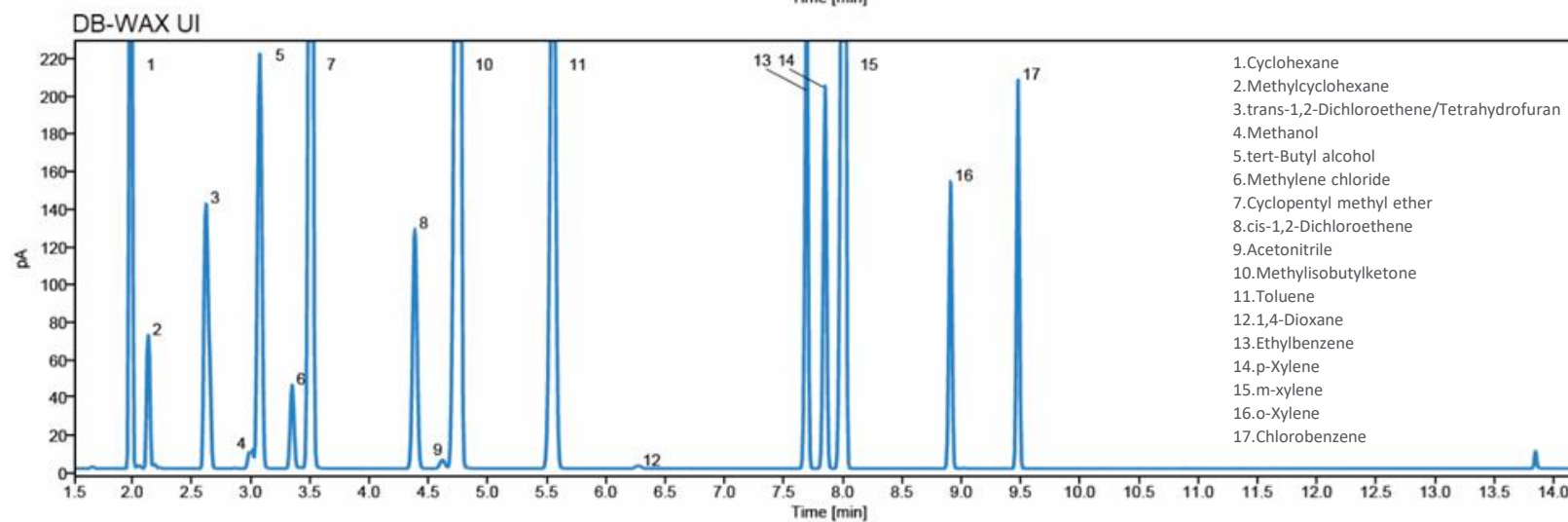


# Class 2A, DB-Select 624 UI & DB-WAX Ultra Inert

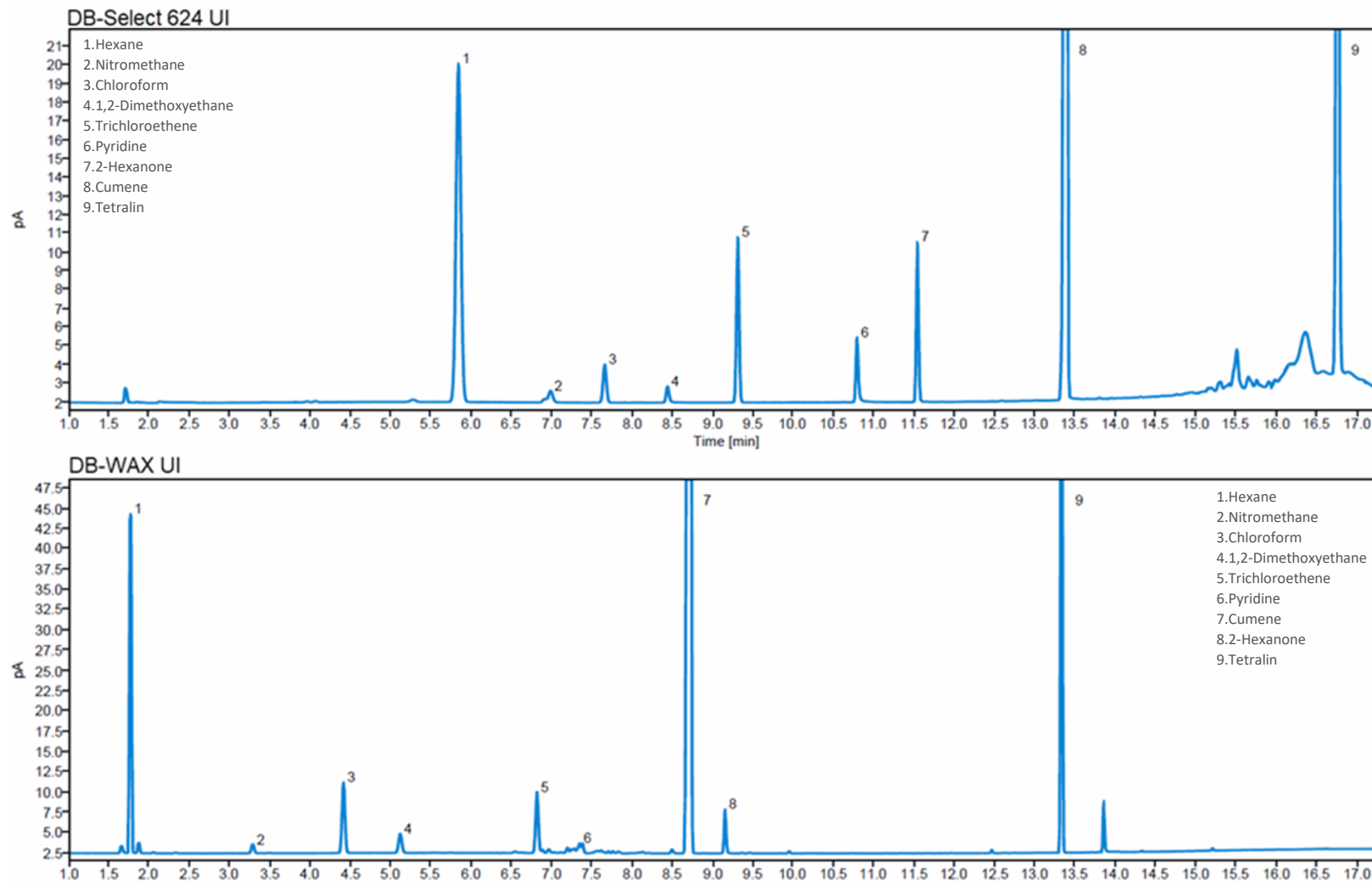
Resolution > 1



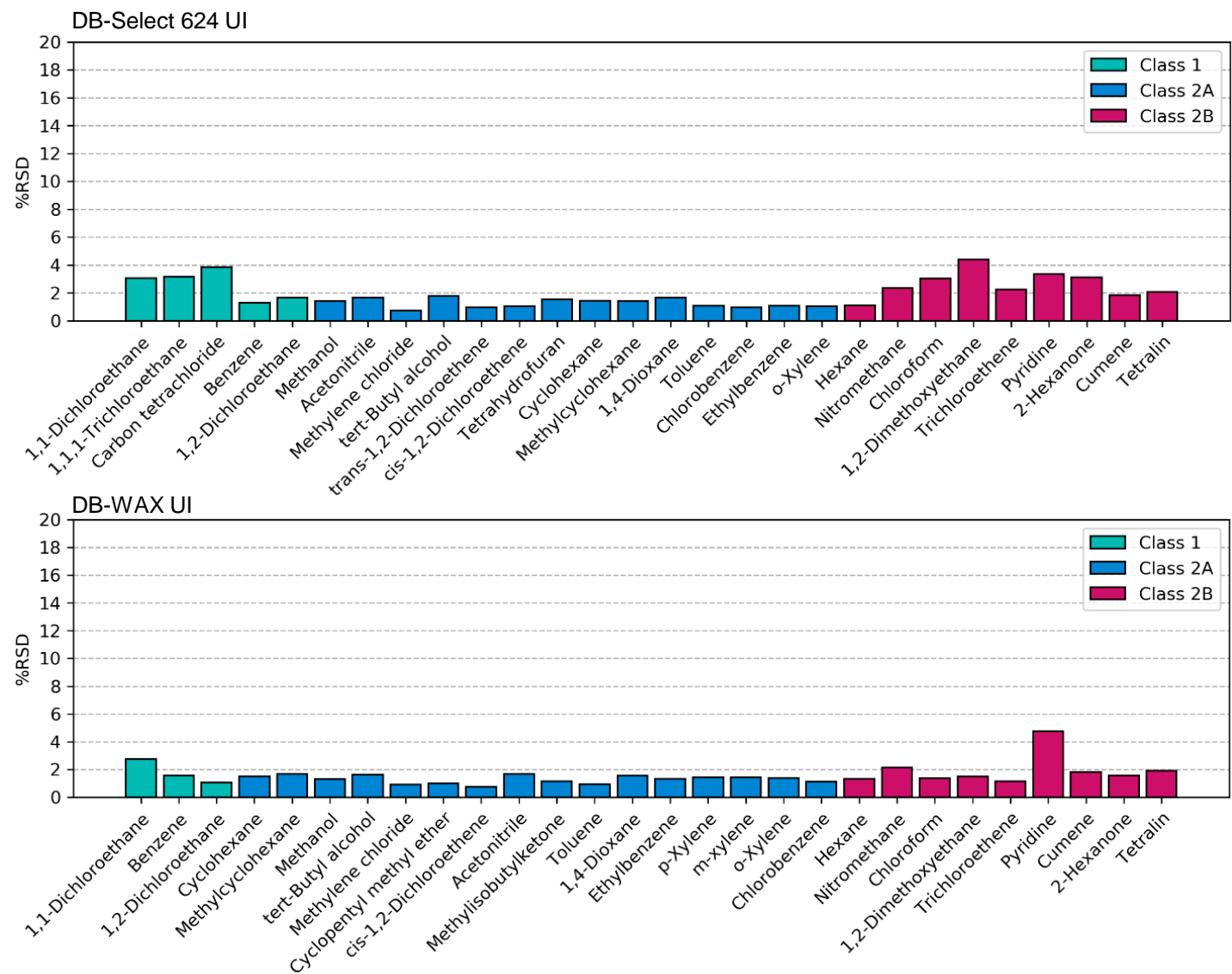
Resolution > 1



# Class 2B, DB-Select 624 UI & DB-WAX Ultra Inert



# Repeatability (n=10)





# USP Method Parameter Comparison for the Agilent 8697 and 7697A Headspace Samplers

Conceptually, both 8897 and 7697A HSSs are very similar, the 8697 HSS can operate with the same method parameters as the 7697A.

Headspace parameters for the USP residual solvents method are shown as an example.

| HS Parameter              | GC/HS            |                  |                  |                  |
|---------------------------|------------------|------------------|------------------|------------------|
|                           | 7890/7697A       | 9000/7697A       | 8890/7697A       | 9000/8697        |
| Sample Loop Volume        | 1 mL             | 1 mL             | 1 mL             | 1 mL             |
| Oven Temperature          | 85 °C            | 85 °C            | 85 °C            | 85 °C            |
| Loop Temperature          | 85 °C            | 85 °C            | 85 °C            | 85 °C            |
| Transfer Line Temperature | 100 °C           | 100 °C           | 100 °C           | 100 °C           |
| Vial Equilibration Time   | 40 min           | 40 min           | 40 min           | 40 min           |
| Injection Duration        | 0.5 min          | 0.5 min          | 0.5 min          | 0.5 min          |
| Vial Size                 | 10 mL            | 10 mL            | 10 mL            | 20 mL*           |
| Vial Shaking              | On, Level 2      | On, Level 2      | On, Level 2      | On, Level 2      |
| Vial Fill Mode            | Flow to pressure | Flow to pressure | Flow to pressure | Flow to pressure |
| Vial Fill Flow            | 50 mL/min        | 50 mL/min        | 50 mL/min        | 50 mL/min        |
| Vial Fill Pressure        | 15 psi           | 15 psi           | 15 psi           | 15 psi           |
| Loop Ramp Rate            | 20 psi/min       | 20 psi/min       | 20 psi/min       | 20 psi/min       |
| Final Loop Pressure       | 0 psi            | 0 psi            | 0 psi            | 4 psi*           |
| Loop Equilibration Time   | 0.05 min         | 0.05 min         | 0.05 min         | 0.05 min         |

\* The most recent application note used 20 mL vials to better accommodate a 6 mL sample volume. Also, the final loop pressure was increased to 4 psi to reduce the potential impact of atmospheric pressure variations, but 0 psi can be used without issue.

# Analysis of Three Classes of Residual Solvents in USP <467> and Chinese Pharmacopoeia by using GC/FID/MSD System



# Configuration Highlights

- This application covers three classes of solvents with a total of up to 62 compounds.
- A purged two-way CFT device was used to split the column effluent 1:1 to the MSD and FID.
- When unknown peaks or unknown solvents appear, this system is the best solution for solvent identification and quantification
- Both MSD and FID signals can be used for quantitative analysis, MSD is a good quantitative supplement for compounds with poor resolution, while FID can expand the linear range.



# Compounds List

The list of compounds in USP <467> and Chinese pharmacopoeia is almost the same.

## Class 2

### Class 1

**Table 1. Class 1 Residual Solvents**  
(Solvents that should be avoided)

| Solvent               | Concentration Limit (ppm) |
|-----------------------|---------------------------|
| Benzene               | 2                         |
| Carbon tetrachloride  | 4                         |
| 1,2-Dichloroethane    | 5                         |
| 1,1-Dichloroethene    | 8                         |
| 1,1,1-Trichloroethane | 1500                      |

### Class 3

**Table 3. Class 3 Residual Solvents**  
(limited by GMP or other quality-based requirements in drug substances, excipients, and drug products)

|                                |                      |
|--------------------------------|----------------------|
| Acetic acid                    | Heptane              |
| Acetone                        | Isobutyl acetate     |
| Anisole                        | Isopropyl acetate    |
| 1-Butanol                      | Methyl acetate       |
| 2-Butanol                      | 3-Methyl-1-butanol   |
| Butyl acetate                  | Methylethylketone    |
| <i>tert</i> -Butylmethyl ether | Methylisobutylketone |
| Cumene                         | 2-Methyl-1-propanol  |
| Dimethyl sulfoxide             | Pentane              |
| Ethanol                        | 1-Pentanol           |
| Ethyl acetate                  | 1-Propanol           |
| Ethyl ether                    | 2-Propanol           |
| Ethyl formate                  | Propyl acetate       |
| Formic acid                    |                      |

Red: liquid injection

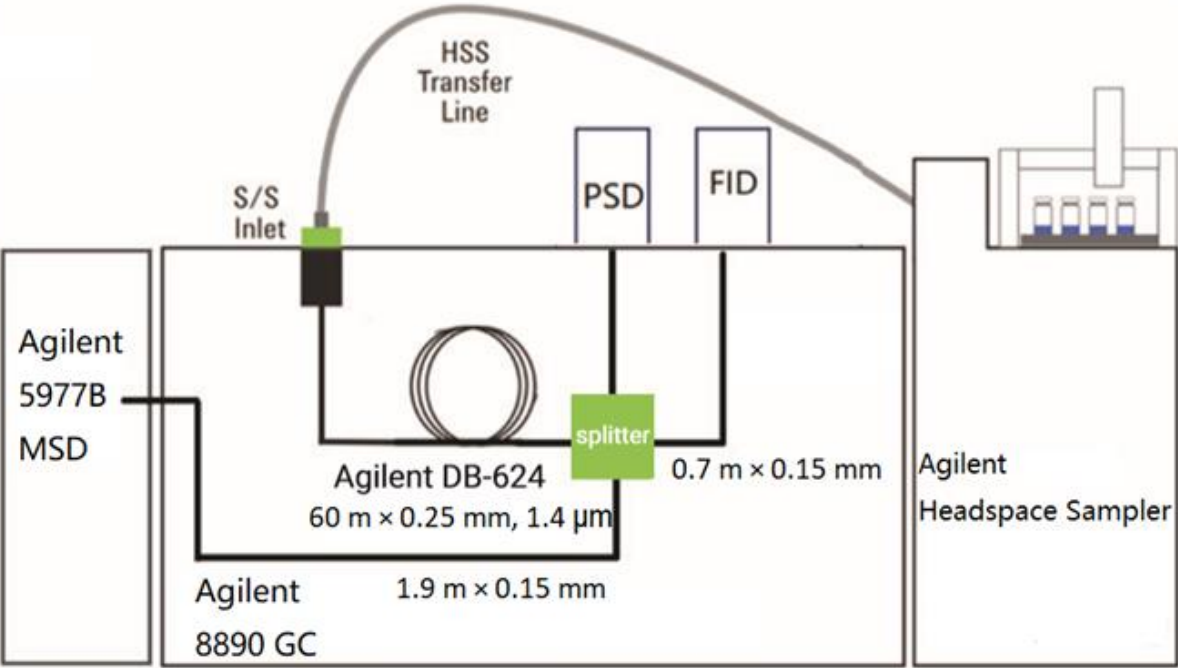
Others: headspace injection

**Table 2. Class 2 Residual Solvents**

| Solvent                       | PDE (mg/day) |
|-------------------------------|--------------|
| Acetonitrile                  | 4.1          |
| Chlorobenzene                 | 3.6          |
| Chloroform                    | 0.6          |
| Cumene                        | 0.7          |
| Cyclohexane                   | 38.8         |
| 1,2-Dichloroethene            | 18.7         |
| 1,2-Dimethoxyethane           | 1.0          |
| <i>N,N</i> -Dimethylacetamide | 10.9         |
| <i>N,N</i> -Dimethylformamide | 8.8          |
| 1,4-Dioxane                   | 3.8          |
| 2-Ethoxyethanol               | 1.6          |
| Ethylene glycol               | 6.2          |
| Formamide                     | 2.2          |
| Hexane                        | 2.9          |
| Methanol                      | 30.0         |
| 2-Methoxyethanol              | 0.5          |
| Methylbutylketone             | 0.5          |
| Methylcyclohexane             | 11.8         |
| Methylene chloride            | 6.0          |
| <i>N</i> -Methylpyrrolidone   | 5.3          |
| Nitromethane                  | 0.5          |
| Pyridine                      | 2.0          |
| Sulfolane                     | 1.6          |
| Tetrahydrofuran               | 7.2          |
| Tetralin                      | 1.0          |
| Toluene                       | 8.9          |
| Trichloroethylene             | 0.8          |
| Xylene*                       | 21.7         |

# Headspace Injection

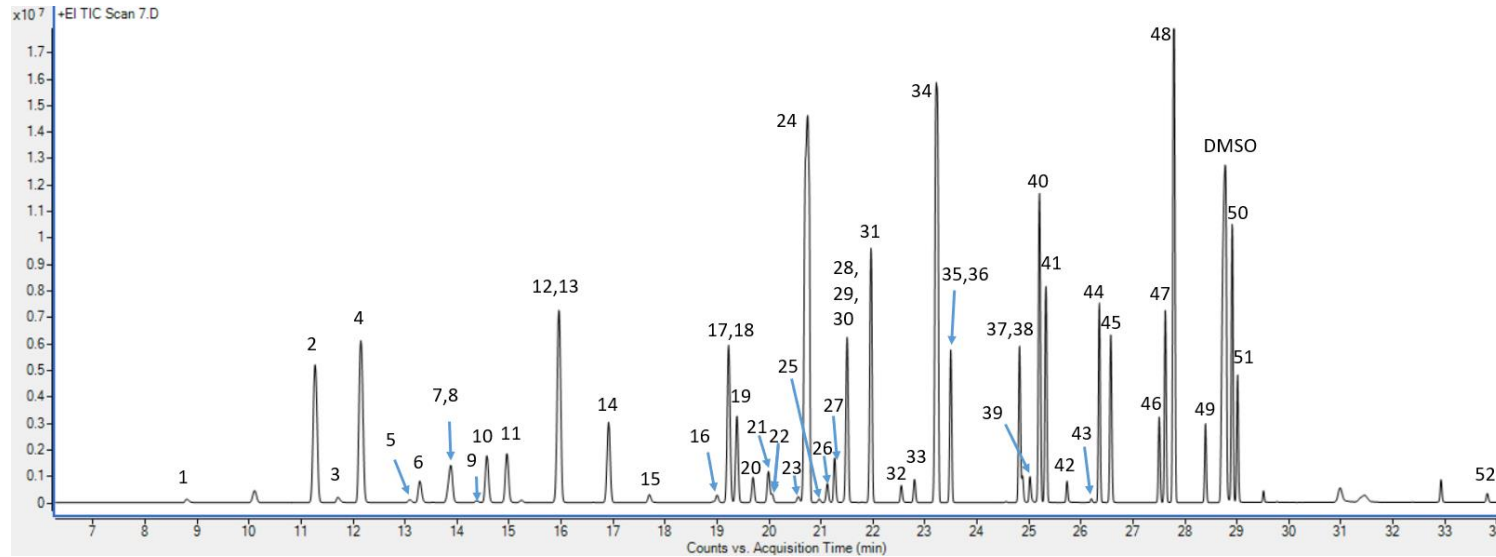
## Instrument conditions



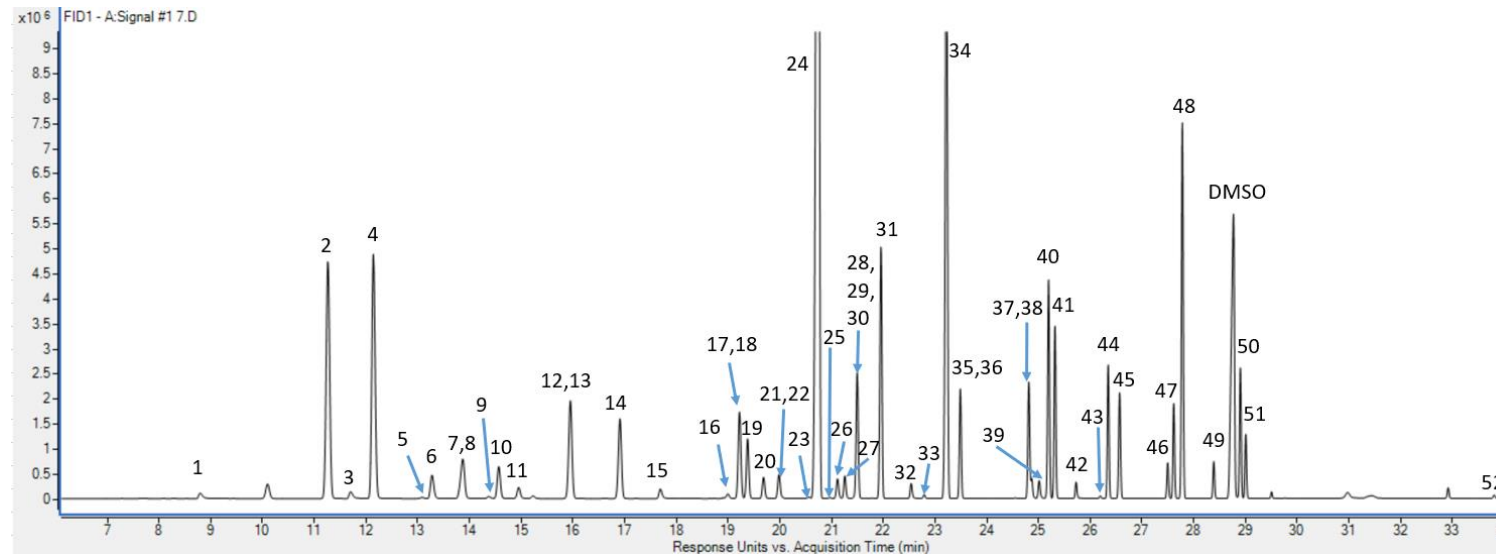
| Agilent 8890 GC           |   |
|---------------------------|---|
| Parameter                 | Value   |
| Inlet                     | SSL, 250 °C, split 10:1   |
| Liner                     | Straight, deactivated, 2 mm ID (part number 5181-8818)                                      |
| CFT Device                | Purged 2-way splitter Split Ratio 1:1 MSD:FID   |
| PSD                       | 3.8 psi constant pressure   |
| Column                    | Agilent DB-624 60 m x 0.25 mm, 1.4 μm (part number 122-1364)                                |
| Carrier                   | Helium, 1 mL/min, constant flow   |
| FID Restrictor            | 0.7 m x 0.15 mm id deactivated fused silica tubing  |
| MSD Restrictor            | 1.9 m x 0.15 mm id deactivated fused silica tubing  |
| Oven                      | 40 °C (10 min), then 5 °C/min to 80 °C, then 12 °C/min to 220 °C (10 min)                   |
| FID                       | Temperature: 250 °C<br>Hydrogen: 30 mL/min<br>Air: 300 mL/min<br>Make-up gas (N2):25 mL/min |
| Transfer line temperature | 250 °C  |
| Agilent 5977B GC/MSD      |   |
| Parameter                 | Value   |
| Ionization type           | EI  |
| Source temperature        | 230 °C  |
| Quad temperature          | 150 °C  |
| Drawout plate             | 3 mm  |
| Tune file                 | Atune.u   |
| Acquisition type          | Scan  |
| Solvent delay             | 6 min   |
| Relative Voltage          | 0   |



# Headspace Injection



MSD scan



FID

# Headspace Injection

## Results of the 52 compounds

| No. | Name                      | RT     | m/z  | Linearity range(µg/mL) | MSD R2 | FID R2 | Area RSD% L4 (n=8) | MDL (MSD) µg/mL L2 (n=8) |
|-----|---------------------------|--------|------|------------------------|--------|--------|--------------------|--------------------------|
| 1   | Methanol                  | 8.818  | 31   | 0.75-150               | 0.9998 | 0.9994 | 2.2                | 0.194                    |
| 2   | Pentane                   | 11.251 | 43   | 0.5-100                | 0.9944 | 0.9997 | 2                  | 0.1428                   |
| 3   | Ethanol                   | 11.73  | 31   | 2-100                  | 0.9999 | 0.9998 | 1.2                | 0.5137                   |
| 4   | Ethyl ether               | 12.142 | 74.1 | 0.5-100                | 0.9911 | 0.9998 | 4.3                | 0.1469                   |
| 5   | 1,1-Dichloroethene        | 13.083 | 61   | 0.004-0.8              | 0.9997 | 0.9986 | 1.7                | 0.0028                   |
| 6   | Acetone                   | 13.283 | 43   | 0.5-100                | 0.9999 | 0.9996 | 2.1                | 0.2265                   |
| 7   | Isopropanol               | 13.854 | 45   | 0.5-100                | 0.9997 | 0.9979 | 2.9                | 0.2446                   |
| 8   | Ethyl formate             | 13.873 | 45   | 0.5-100                |        |        | 4.3                | 0.2449                   |
| 9   | Acetonitrile              | 14.39  | 41   | 0.1-20                 | 0.9996 | 0.9984 | 4.2                | 0.0319                   |
| 10  | Methyl acetate            | 14.564 | 43   | 0.5-100                | 0.9998 | 0.9998 | 2.7                | 0.4236                   |
| 11  | Methylene chloride        | 14.947 | 84   | 0.15-30                | 0.9997 | 0.9997 | 2.1                | 0.0326                   |
| 12  | 2-Methoxy-2-methylpropane | 15.938 | 73   | 0.1-20                 | 0.9988 | 0.9998 | 2.1                | 0.0352                   |
| 13  | trans-1,2-Dichloroethene  | 15.979 | 95.9 | 0.235-47               | 0.9969 | 0.9998 | 1.7                | 0.065                    |
| 14  | Hexane                    | 16.899 | 57   | 0.1-20                 | 0.9995 | 0.9998 | 2.2                | 0.0739                   |
| 15  | 1-Propanol                | 17.712 | 31   | 0.5-100                | 0.9995 | 0.9996 | 2                  | 0.1799                   |
| 16  | Nitromethane              | 19     | 46   | 0.5-100                | 0.9999 | 0.9991 | 1.9                | 0.2521                   |
| 17  | cis-1,2-Dichloroethene    | 19.21  | 96   | 0.235-47               | 0.9988 | 0.9999 | 2.5                | 0.0447                   |
| 18  | 2-Butanone                | 19.225 | 43   | 0.5-100                | 0.998  | 0.9999 | 2.3                | 0.1471                   |
| 19  | Ethyl acetate             | 19.375 | 43   | 0.5-100                | 0.9986 | 0.9997 | 1.4                | 0.3054                   |
| 20  | 2-Butanol                 | 19.688 | 45   | 0.5-100                | 0.9998 | 0.9999 | 2.4                | 0.2371                   |
| 21  | Tetrahydrofuran           | 19.985 | 42   | 0.18-36                | 0.9998 | 0.9998 | 2.1                | 0.0532                   |
| 22  | Chloroform                | 20.054 | 83   | 0.015-3                | 0.9997 |        | 1.6                | 0.0058                   |
| 23  | 1,1,1-Trichloroethane     | 20.546 | 97   | 0.005-1                | 0.9999 | 0.9998 | 1.3                | 0.0025                   |
| 24  | Cyclohexane               | 20.707 | 84   | 1.0-49 (195)*          | 0.9908 | 0.9997 | 1.8                | 0.188                    |
| 25  | Carbon tetrachloride      | 20.962 | 117  | 0.002-0.4              | 0.9998 | 0.9992 | 2.8                | 0.002                    |

| No. | Name                          | RT     | m/z  | Linearity range(µg/mL) | MSD R2 | FID R2 | Area RSD% L4 (n=8) | MDL (MSD) µg/mL L2 (n=8) |
|-----|-------------------------------|--------|------|------------------------|--------|--------|--------------------|--------------------------|
| 27  | 1,2-Dimethoxyethane           | 21.265 | 45   | 0.5-100                | 0.9999 | 0.9995 | 1                  | 0.2561                   |
| 28  | Benzene                       | 21.442 | 78   | 0.001-0.2              | 0.9995 | 0.9998 | 5.8                | 0.0008                   |
| 29  | 1,2-Dichloroethane            | 21.442 | 62   | 0.01-0.5               | 0.9989 |        | 1.5                | 0.0016                   |
| 30  | Isopropyl acetate             | 21.496 | 61   | 0.5-100                | 0.9985 | 0.9998 | 0.8                | 0.1636                   |
| 31  | Heptane                       | 21.956 | 71   | 0.1-20                 | 0.9974 | 0.9996 | 2.4                | 0.0343                   |
| 32  | 1-Butanol                     | 22.547 | 56   | 0.5-100                | 0.9994 | 0.9998 | 2.4                | 0.1717                   |
| 33  | Trichloroethylene             | 22.791 | 130  | 0.015-3                | 0.9999 | 0.9999 | 1.8                | 0.0065                   |
| 34  | Methyl cyclohexane            | 23.208 | 83   | 0.3-15 (59)*           | 0.9989 | 0.9997 | 2.3                | 0.0722                   |
| 35  | 1,4-Dioxane                   | 23.489 | 88   | 0.095-19               | 0.9999 | 0.9999 | 3.3                | 0.0549                   |
| 36  | Propyl acetate                | 23.491 | 43   | 0.5-100                | 0.9966 |        | 3                  | 0.2675                   |
| 37  | 4-Methyl-2-pentanone          | 24.815 | 43   | 0.5-100                | 0.9985 | 0.9999 | 2.2                | 0.1429                   |
| 38  | Isoamyl alcohol               | 24.879 | 55.1 | 0.5-100                | 0.9991 | 0.9996 | 2.4                | 0.2562                   |
| 39  | Pyridine                      | 25.024 | 79   | 2-100                  | 0.9992 | 0.9997 | 2.1                | 0.5016                   |
| 40  | Toluene                       | 25.196 | 91   | 0.225-22 (44)*         | 0.9964 | 0.9998 | 2.1                | 0.0651                   |
| 41  | Isobutyl acetate              | 25.322 | 56   | 0.5-100                | 0.9958 | 0.9999 | 2.1                | 0.1784                   |
| 42  | 1-Pentanol                    | 25.735 | 42   | 0.5-100                | 0.9996 | 0.9998 | 2.1                | 0.3319                   |
| 43  | 2-Hexanone                    | 26.201 | 58   | 0.06-3                 | 0.9995 | 0.9998 | 2.1                | 0.0107                   |
| 44  | Butyl acetate                 | 26.351 | 43   | 0.5-100                | 0.9957 | 0.9999 | 2.3                | 0.2502                   |
| 45  | Tetrahydrothiophene           | 26.571 | 88   | 0.5-100                | 0.9996 | 0.9999 | 1.4                | 0.18                     |
| 46  | Chlorobenzene                 | 27.503 | 112  | 0.09-18                | 0.9999 | 0.9997 | 2.5                | 0.0215                   |
| 47  | Ethylbenzene                  | 27.618 | 91   | 0.09-18                | 0.9986 | 0.9997 | 4.1                | 0.0288                   |
| 48  | m,p-xylene                    | 27.782 | 106  | 0.4-40 (80)*           | 0.9963 | 0.9997 | 3.3                | 0.1074                   |
| 49  | o-xylene                      | 28.393 | 91   | 0.05-10                | 0.9999 | 0.9996 | 2.6                | 0.0173                   |
| 50  | Isopropylbenzene              | 28.904 | 105  | 0.1-20                 | 0.9983 | 0.9996 | 2.4                | 0.0391                   |
| 51  | Anisole                       | 29.011 | 108  | 0.5-100                | 0.9999 | 0.9997 | 2.8                | 0.1892                   |
| 52  | 1,2,3,4-Tetrahydronaphthalene | 33.814 | 104  | 0.015-3                | 0.9998 | 0.9993 | 2                  | 0.0045                   |

# Liquid Injection

## Compounds list in liquid injection

2-Methoxyethanol

2-Ethoxyethanol

N,N-dimethylformamide

N,N-dimethylacetamide

Acetic acid

Formic acid

Ethylene glycol

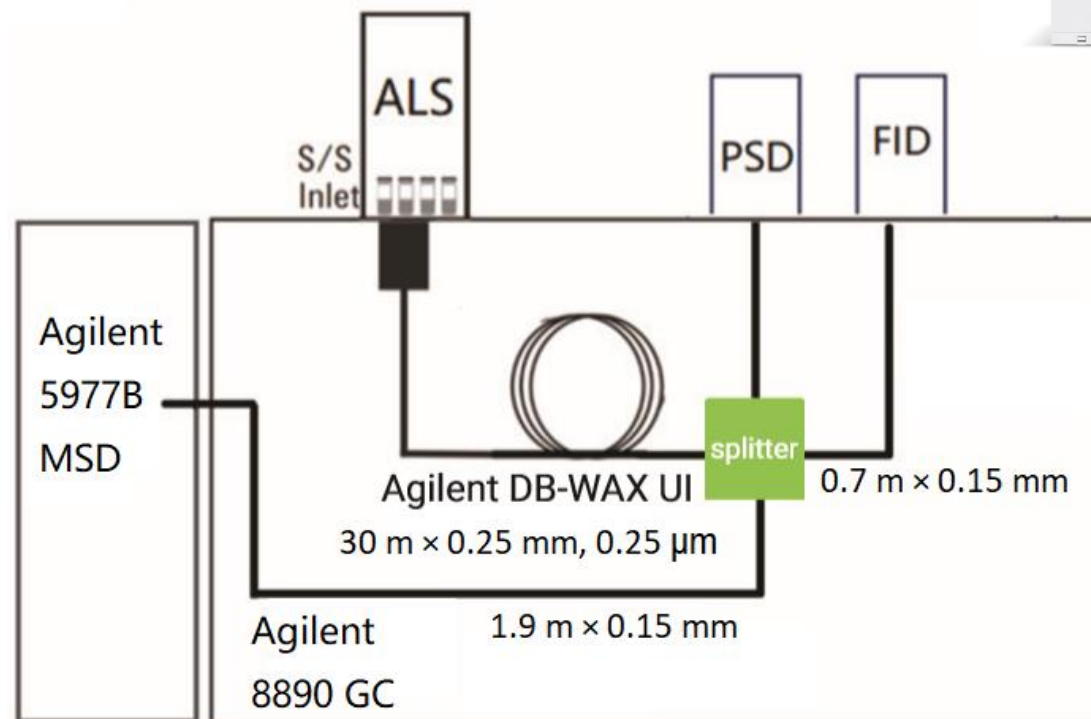
N-methylpyrrolidone

Formamide

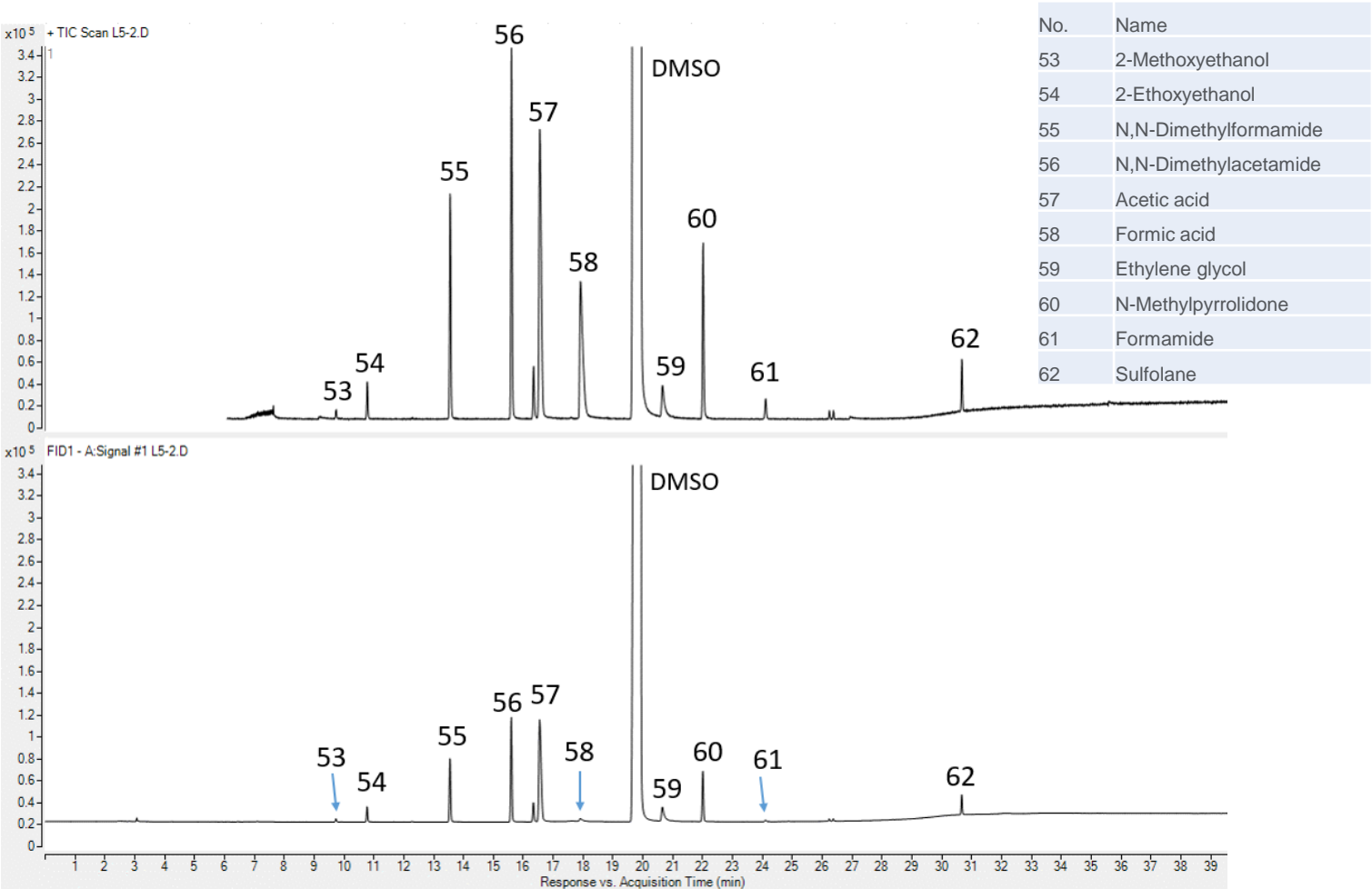
Sulfolane

Standards: USP 467 Class 2C (Agilent p/n: 5190-0493)  
acetic acid (99.8%, purity), formic acid (98%, purity)

## Instrument configuration



# Liquid Injection



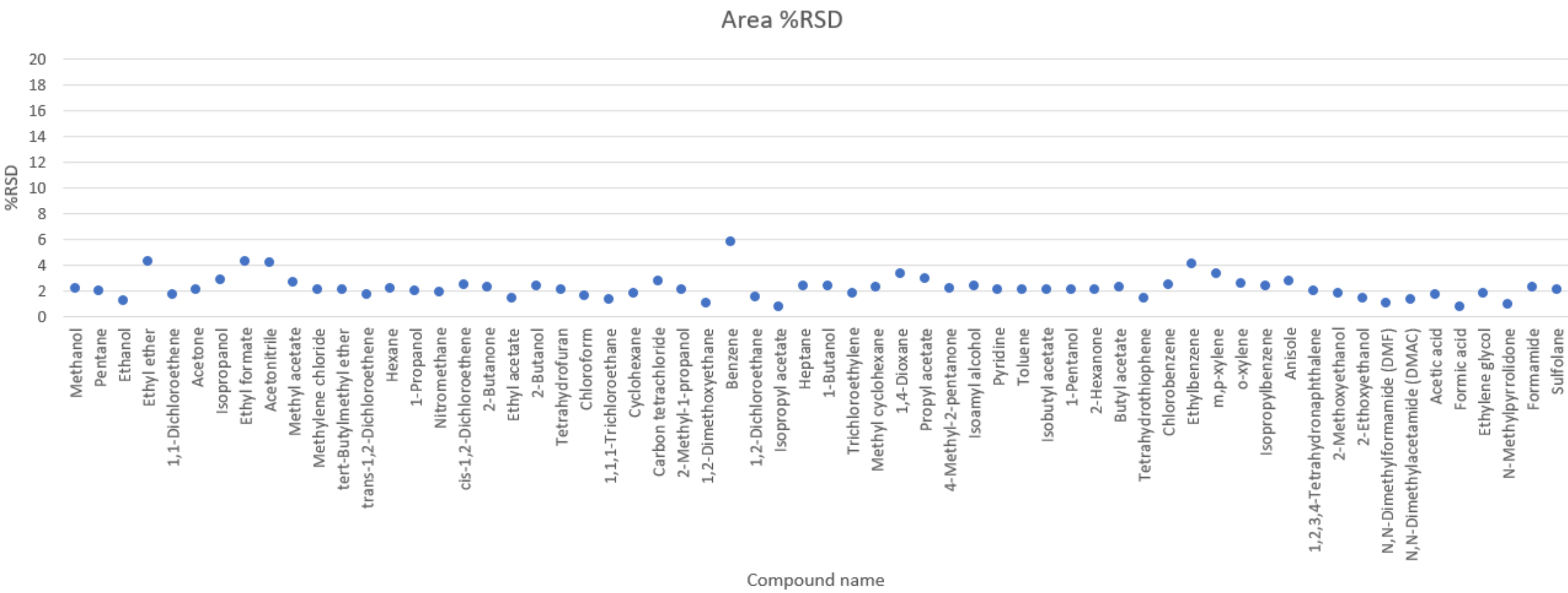
| Agilent 8890 GC           |  |
|---------------------------|--|
| Parameter                 | Value  |
| Inlet                     | SSL, 250 °C, split 30:1  |
| Liner                     | Ultra Inert, split, low pressure drop, glass wool (p/n: 5190-2295)                           |
| Injection volume          | 0.5 uL   |
| CFT Device                | Purged 2-way splitter Split Ratio 1:1 MSD:FID  |
| PSD                       | 3.8 psi constant pressure  |
| Column                    | Agilent DB-wax UI 30 m x 0.25 mm, 0.25 µm (part number 122-7032UI)                           |
| Carrier                   | Helium, 1 mL/min, constant flow  |
| FID Restrictor            | 0.7 m x 0.15 mm id deactivated fused silica tubing   |
| MSD Restrictor            | 1.9 m x 0.15 mm id deactivated fused silica tubing   |
| Oven                      | 40 °C, then 5 °C/min to 160 °C, then 10 °C/min to 220 °C (10 min)                            |
| FID                       | Temperature: 250 °C<br>Hydrogen: 30 mL/min<br>Air: 300 mL/min<br>Make-up gas (N2): 25 mL/min |
| Transfer line temperature | 250 °C   |
| Agilent 5977B GC/MSD      |  |
| Parameter                 | Value  |
| Ionization type           | EI   |
| Source temperature        | 230 °C   |
| Quad temperature          | 150 °C   |
| Drawout plate             | 3 mm   |
| Tune file                 | Atune.u  |
| Acquisition type          | Scan   |
| Solvent delay             | 6 min  |
| Relative Voltage          | 0  |

# Liquid Injection

## Results of the 10 compounds

| No. | Name                         | RT     | m/z | Linearity range<br>μg/mL | R <sup>2</sup> |        | Area RSD%<br>L4 (n=8) | MDL<br>(MSD)<br>μg/mL |
|-----|------------------------------|--------|-----|--------------------------|----------------|--------|-----------------------|-----------------------|
|     |                              |        |     |                          | MSD            | FID    |                       |                       |
| 53  | 2-Methoxyethanol             | 9.783  | 45  | 5-50                     | 0.9984         | 0.9995 | 1.8                   | 0.68                  |
| 54  | 2-Ethoxyethanol              | 10.816 | 59  | 16-161                   | 0.9973         | 0.9987 | 1.4                   | 1.93                  |
| 55  | N,N-Dimethylformamide (DMF)  | 13.607 | 73  | 88.3-883                 | 0.9997         | 0.9999 | 1                     | 2.19                  |
| 56  | N,N-Dimethylacetamide (DMAC) | 15.667 | 87  | 109.4-1094               | 0.9997         | 0.9996 | 1.3                   | 2.58                  |
| 57  | Acetic acid                  | 16.493 | 60  | 400-3000                 | 0.9984         | 0.9997 | 1.7                   | 90.12                 |
| 58  | Formic acid                  | 17.774 | 46  | 400-3000                 | 0.9995         | 0.9939 | 0.8                   | 120                   |
| 59  | Ethylene glycol              | 20.652 | 31  | 62.2-622                 | 0.9983         | 0.9982 | 1.8                   | 4.44                  |
| 60  | N-Methylpyrrolidone          | 22.074 | 98  | 53-530                   | 0.9995         | 0.9997 | 0.9                   | 3.02                  |
| 61  | Formamide                    | 24.157 | 45  | 22-221                   | 0.9992         | 0.9986 | 2.3                   | 2.11                  |
| 62  | Sulfolane                    | 30.706 | 120 | 16-160                   | 0.9994         | 0.9997 | 2.1                   | 1.33                  |

# Repeatability (n=8) for 62 Compounds





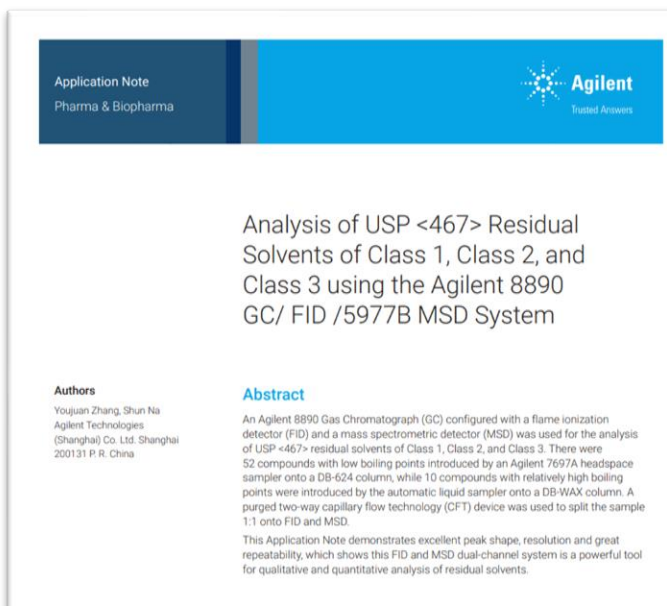
# Summary

- Residual solvents of Class 1, 2, and 3 were tested using the Agilent 8890 GC/FID/MSD system.
- For new drug development and quality control, FID and MSD dual-channel configurations can be powerful tools for solvent residue analysis.
- MSD analysis can avoid the uncertainty of more than 60 solvents involved in drug production.
- When unknown peaks or unknown solvents appear, this system is the best solution for solvent identification and quantification

## Application Notes

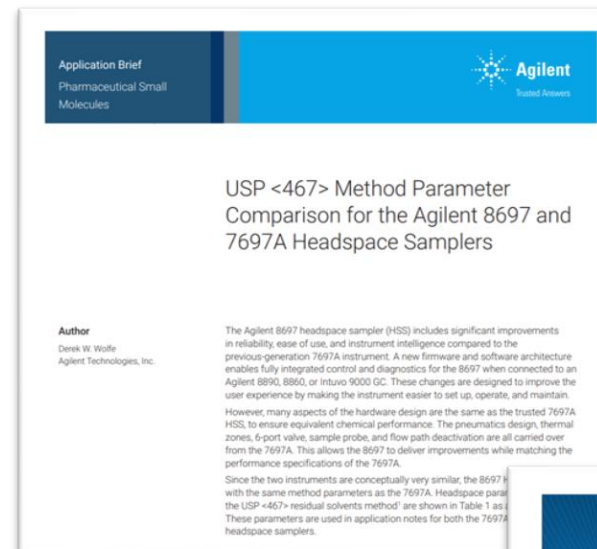


[Residual Solvents Analysis Using an Agilent Intuvo 9000 GC with 8697 Headspace Sampler](#)



[Analysis of USP <467> Residual Solvents of Class 1, Class 2, and Class 3 using the Agilent 8890 GC/ FID /5977B MSD System](#)

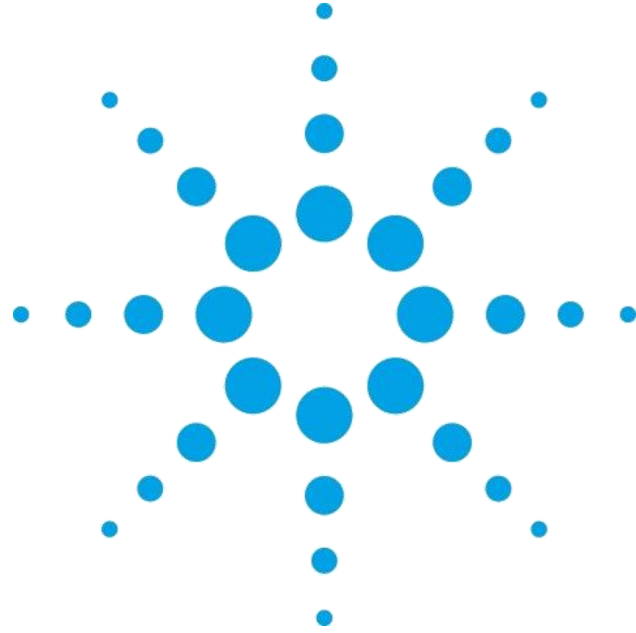
## Application Brief



[USP <467> Method Parameter Comparison for the Agilent 8697 and 7697A Headspace Samplers](#)



<https://www.agilent.com/cs/library/application/s/5991-8032EN.pdf>



# Thank you for your attention!