

Analysis of Extractables & Leachables in Pharmaceutical Products



Regulatory & Analytical Aspects

Dr. Andreas Tei
Global Pharma Segment Manager
Agilent Technologies

Outline

Section #1

- Introduction
- Defining extractables and leachables
- Guidelines
- Effects on biologic drug products
- The four essential steps of a study

Section #2

- Application Example: analysis of elemental E&Ls by ICP-MS
- Application Example: analysis of volatile E&Ls by GC-MS
- Application Example: analysis of non volatile E&Ls by LC-MS
- Appendix / References

Introduction

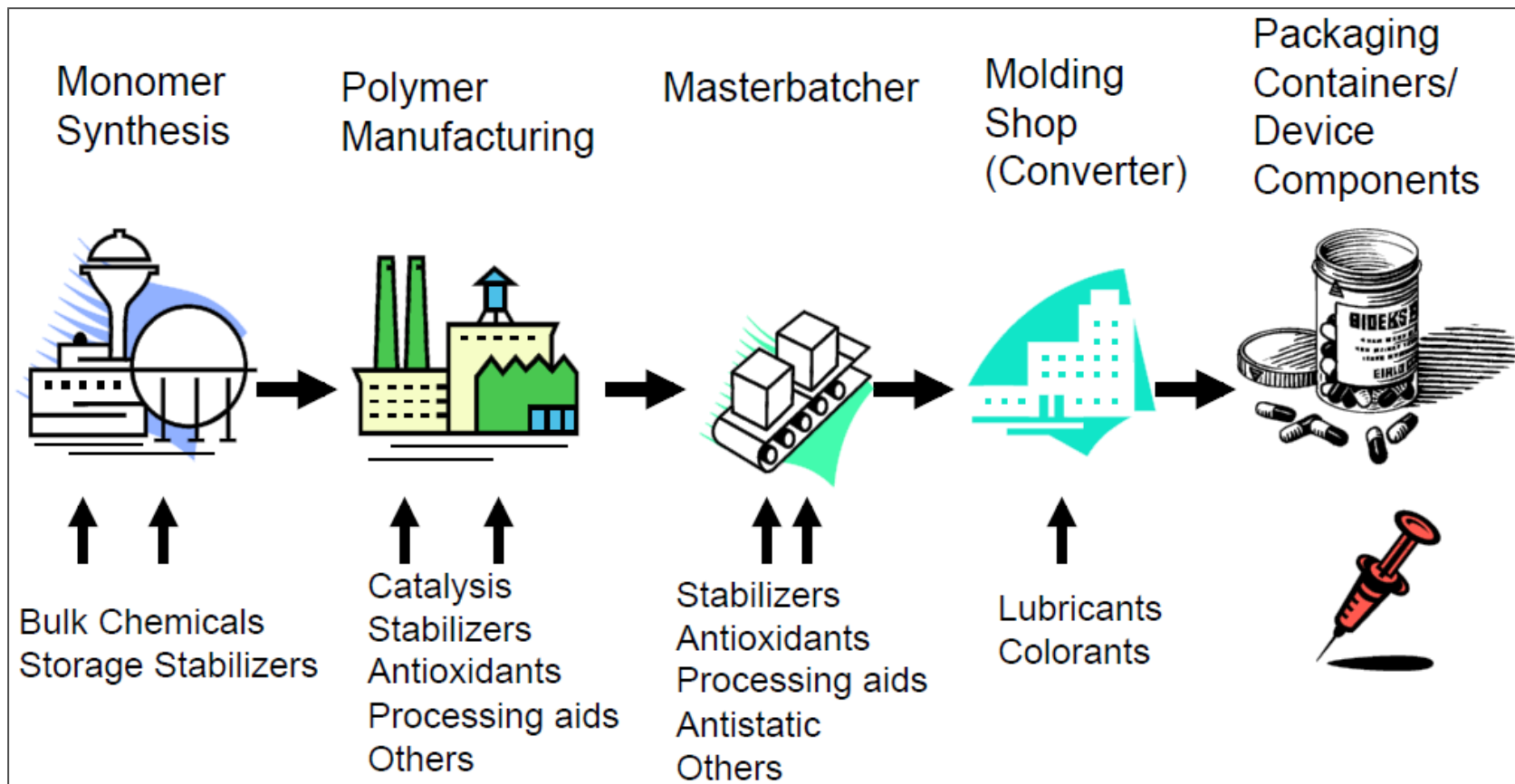
Drug Containers And Modern Drug Delivery Systems

Drug containers and modern drug delivery systems meant to protect a drug from environmental contamination but they are actually themselves a source of contamination



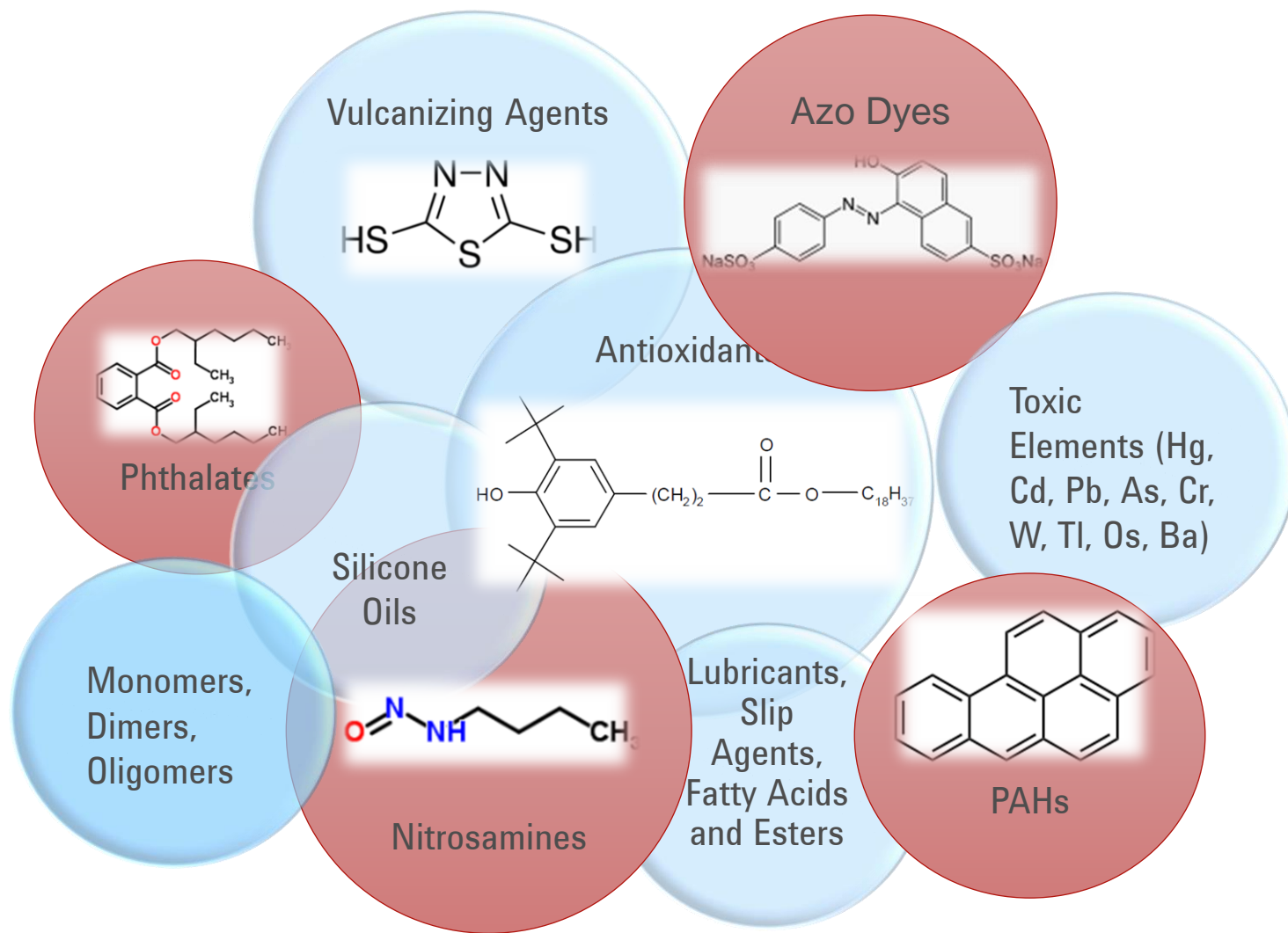
Plastic Materials: Source of Contamination

Sources of extractables are plastic and elastomeric components (monomers, polymeric initiators, plasticizers, etc.) ink and adhesives (label) and degradation products (processing, storage, sterilization)



Cindy Zweiben, Pfizer, Inc., Characterization of Extractables and Leachable in Parenteral Drug Products

Compounds Identified as Extractables/Leachables

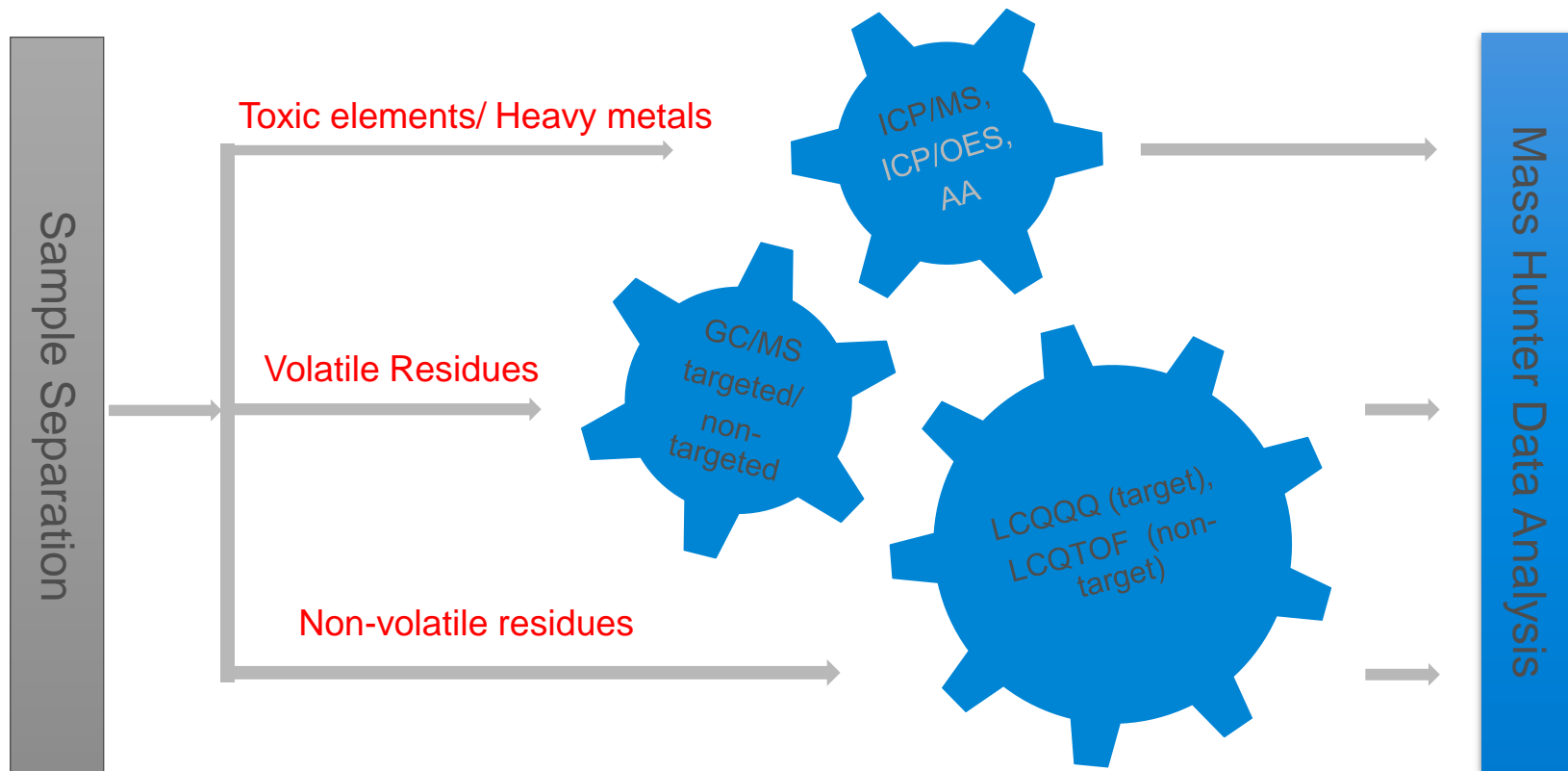


Wide variety of Chemical Classes, Polarity, Molecular Weights, Properties

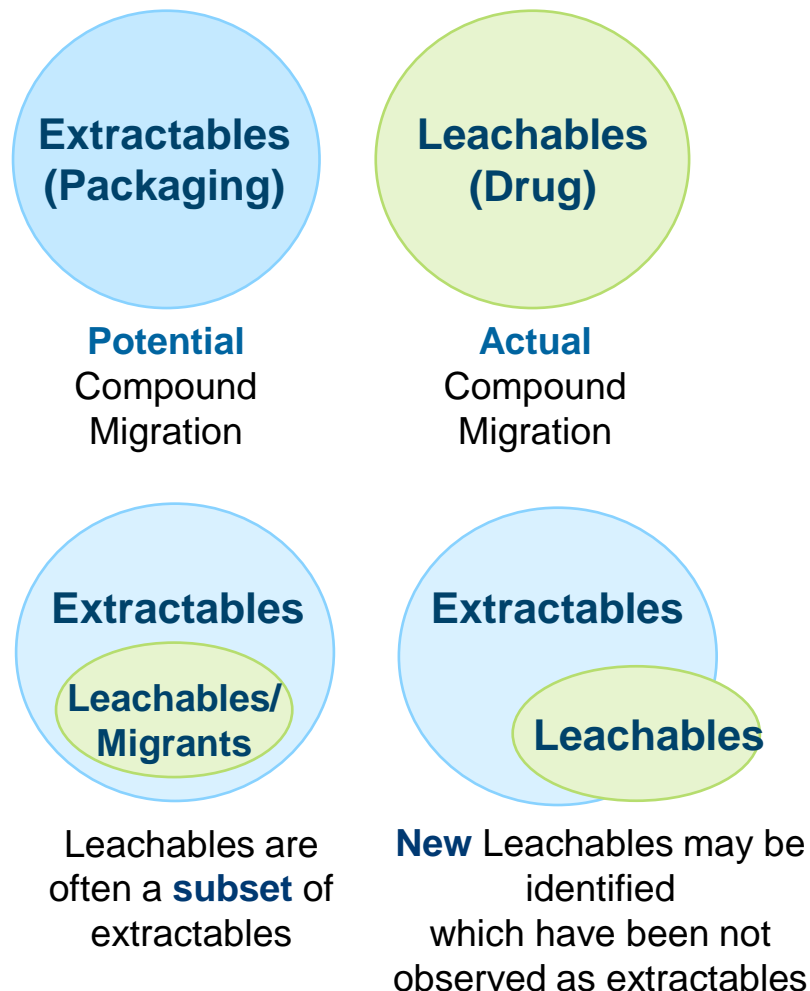
Various Analytical Technologies Required

Agilent delivers the most comprehensive analytical solutions portfolio

Objective: To detect a wide class of known and unknown organic/inorganic compounds that maybe present in container closure systems at levels links to risk assessment threshold levels



Defining Extractables, Leachables, Migrants



Extractable

Chemical compounds that can be extracted out of packaging component

- Analyze **packaging component** at
 - High-temperatures**: to obtain the worst case leachable profile
 - Solvent extraction**: polar and non-polar solvent to mimic similar properties as drug product

Leachable

- Chemical compounds from packaging component that leach into the drug product
- Analyze **drug product** at
 - Normal conditions
 - Simulate extended storage conditions

Migrants

- Crossed the primary packaging material barrier from secondary and tertiary packaging, accumulating in the drug product



FDA Regulation For Container Closure Systems (CCS)

U.S. FDA 21 CFR 211.94(a) statement (April 2015)

“(a) Drug product containers and closures shall not be reactive, additive, or absorptive so as to alter the safety, identity, strength, quality, or purity of the drug beyond the official or established requirements.

...(d) Standards or specifications, methods of testing, and, where indicated, methods of cleaning, sterilizing, and processing to remove pyrogenic properties shall be written and followed for drug product containers and closures.”

<http://www.accessdata.fda.gov/scripts/cdrh/cfdocs/cfcfr/CFRSearch.cfm?fr=211.94>



Why Worry about Extractables/Leachables ?

2014: FDA Drug Recalls Surges over 836 in 2014!

2014: FDA data shows the last two years have seen almost as many recalls (2,061) as the previous nine years combined (2,217)—and that's only counting the first seven months of 2014. *Ref: raps.org August 2014*

August 2015: FDA warns against use of Becton-Dickinson (BD) 3 ml and 5 ml

Syringes: <http://www.fda.gov/Safety/MedWatch/SafetyInformation/SafetyAlertsforHumanMedicalProducts/ucm458955.htm>

Sodium Valproate Zentiva 500mg Gastro-resistant Tablets, PL 17780/0454, Zentiva Livery

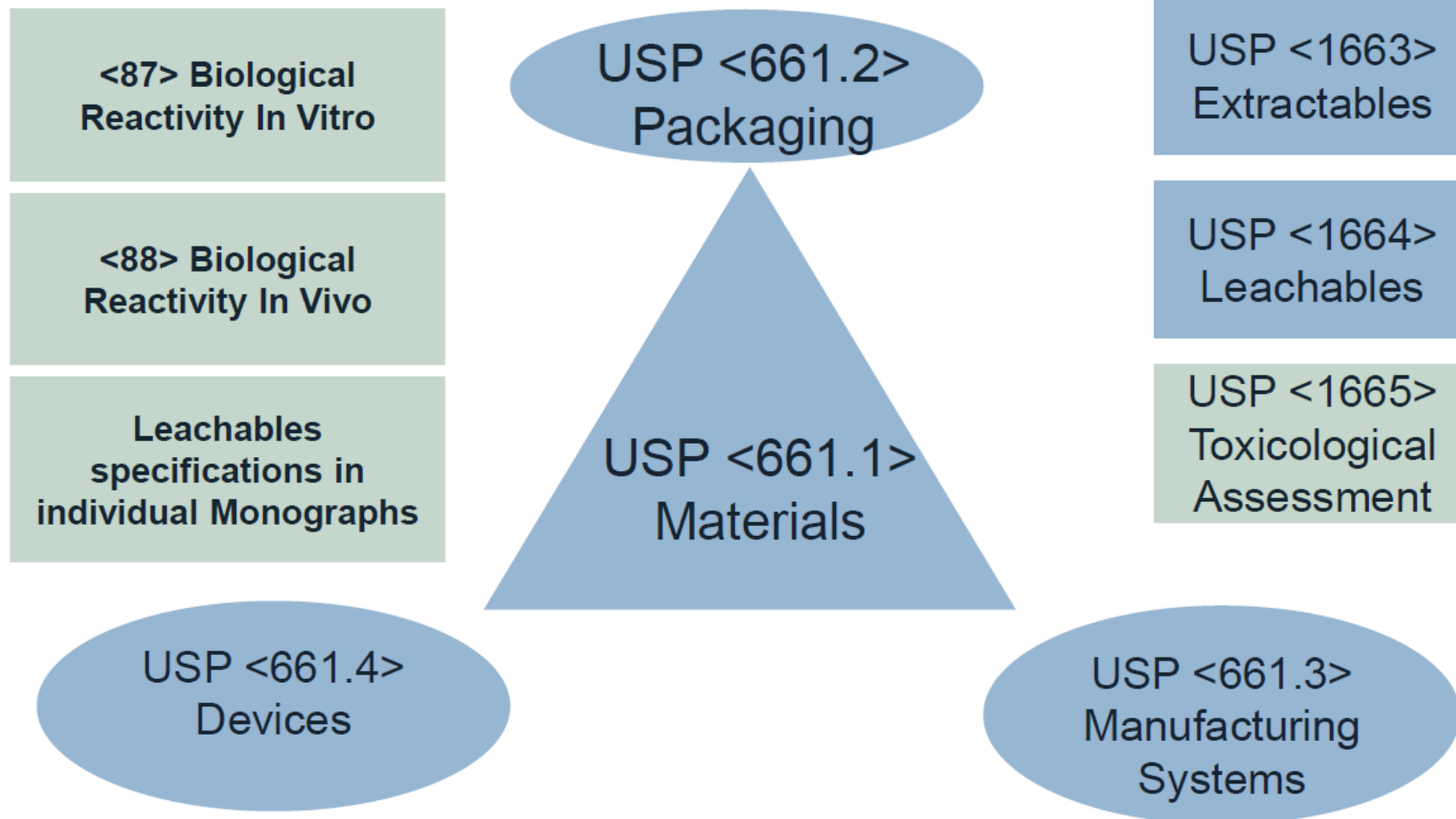
Batch Number	Expiry date	Pack Size	First Distributed
J601	Aug 2016	1 x 100	01 Oct 2014

We have been made aware of an unusual odour associated with the above batches of Epilim / Sodium Valproate Gastro-resistant tablets. The odour has variously been described as 'fishy', 'sweaty armpits' and 'cannabis'. An investigation by Sanofi has identified the root cause as the aluminium foil used in the packaging of affected batches. No impact on the quality or efficacy of the tablets has been detected. Sanofi has now ceased using this foil supplier.

https://assets.digital.cabinet-office.gov.uk/media/55191d6fe5274a142e000069/EL__15_A_02.pdf



USP Chapters dealing with E&L



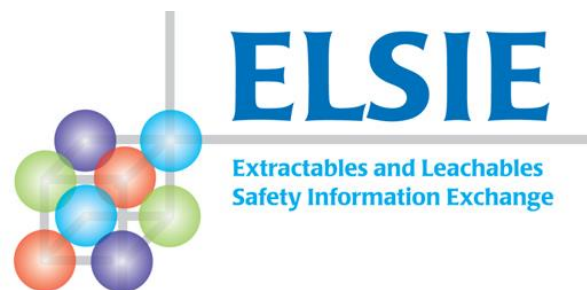
Ref: Denise R. Jenke, Daniel L. Norwood, and Desmond G Hunt



Guidelines Delivered By Pharma Industry Expert Working Groups



PQRI (Product Quality Research Institute) is a working group established to develop **regulatory guidance** for Extractable/Leachable analysis, which is also recognized by the FDA



PQRI guidance for OINDP

(Orally inhaled and nasal drug products)

PQRI issued guidance for OINDP:

“Safety thresholds and best practices for extractables and leachables in orally inhaled and nasal drug products (OINDP) also Applicable to parenteral and injectable products (PODP)

PQRI established safety thresholds for leachables:

- **Safety Concern Threshold (SCT)** $\leq 0.15 \mu\text{g/day}$ patient exposure which species represent **no risk**
- **Qualification Threshold (QT)** $\leq 5 \mu\text{g/day}$ patient exposure which a leachable is **not considered for safety qualification**. **Lower** threshold applies to PAH's, nitrosamines, and 2-mercaptobenzothiazole
- **Estimated Analytical Evaluation Threshold (AET) ($\mu\text{g/g}$)** = $(\text{SCT} \times \text{total labeled doses}) / (\text{Doses per day} \times \text{mass of component})$

“Best practices” include controlled extraction studies and leachables studies.



Inorganic Impurities

New USP general chapters <232> and <233> for elemental impurities



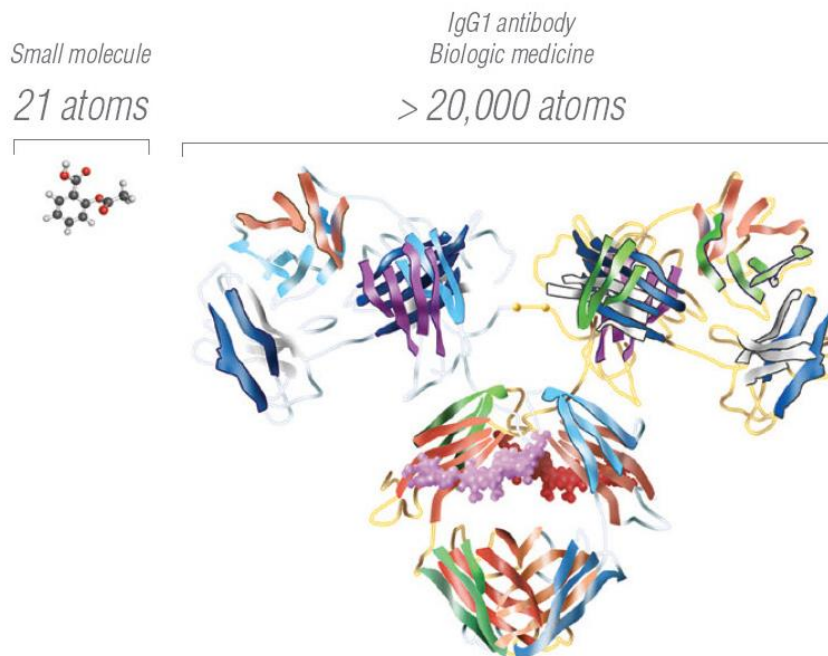
- Reagents, Ligands, Catalysts
- Manufacturing Aids
- Inks and Dyes

USP<232> defines the analyte limits, while **USP<233>** defines sample preparation options including closed vessel microwave digestion, and recommends the use of modern instrumentation, such as multi-element ICP-MS and ICP-OES techniques. Analytical equipment qualification under USP<233> is based on performance testing, and includes requirements to demonstrate accuracy, repeatability, and the unequivocal identification of analytes.

Special Concerns About E&L Effects on Biologics

Even contaminations at trace levels with reactive E&L impurities can be deleterious for protein based drugs and will cause severe harm for the patient's health (immunogenic reaction)

E&L compound
as Impurity



Contact materials are: plastics/elastomers, glass and stainless steel surfaces



Sources of E&Ls as contaminants in biological drugs

☐ Plastics / Elastomers

- Crosslinking agents, volatile organic E&Ls

☐ Glass surfaces

- Al^{3+} ; $\text{Fe}^{3+/2+}$; Ca^{2+} ; Ba^{2+} ; Mn^{2+} ; Zn^{2+}

☐ Stainless steel surfaces

- $\text{W}^{6+/4+}$; $\text{Fe}^{3+/2+}$; $\text{Cr}^{3+/2+}$; Ni^{2+}

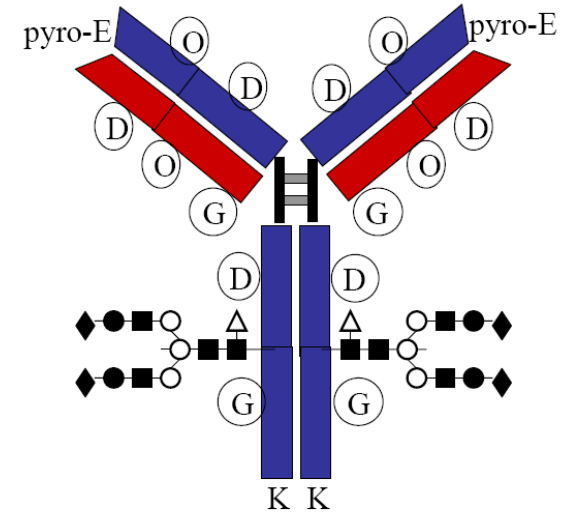
☐ Sterile filtration processes are often a source of contamination

☐ Sterilization processes (steam autoclaving / gamma radiation) of drug containers will affect the concentration of leachables within the drug product

☐ Complexing agents (EDTA) facilitate migration of metal ions

E&Ls may affect protein drug products by:

- Aggregation
- Increase in particulates*
- Oxidation
- Unfolding
- Formation of clipped variants
- Formation of Protein Adducts
- Post translational events during fermentation (glycosylation)
- Altered protein translation

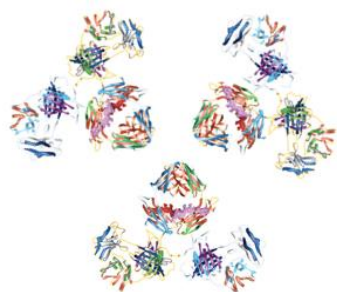


* See also new USP monograph <787> PARTICULATE MATTER IN THERAPEUTIC INJECTIONS

Ref: Ingrid Markovic, CBER Presentation USP/PQRI E/L Workshop April 2014
https://www.usp.org/sites/default/files/usp_pdf/EN/meetings/09_markovich_presentation.pdf

Critical Quality Attributes & Testing Methods for mAbs

Aggregate Analysis of Monoclonal Antibody



SEC
Aggregation

Disulfide Shuffling **RP**

Pyro-Glutamate **IEC**

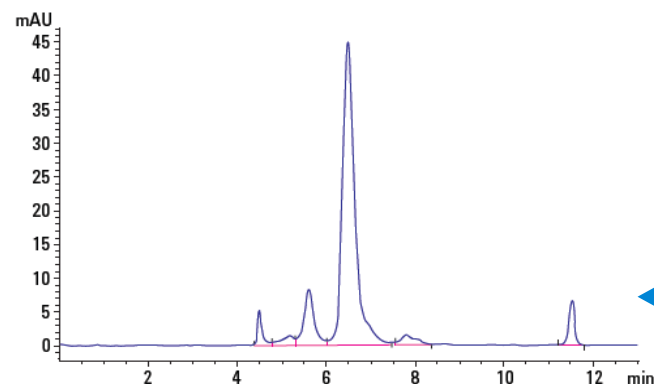
Deamidation /Oxidation **IEC**

Fragmentation (Hinge) **RP**

Truncation (Lys 0, 1, 2) **IEC**

Glycosylation (G0, G1, G2) **HILIC**

Column: Agilent AdvanceBio SEC 300Å, 2.7 µm, 7.8 x (p/n PL1180-5301)
Flow rate: 1 mL/min
Mobile phase: 150 mM phosphate buffer, pH 7.0
Wavelength: 220 nm
Temperature: ambient
Injection volume: 5 µL
Sample: IgG



Agilent Technologies

Key Principles of an E&L Study

1

Evaluating The Interactions Between Packaging Material And The Pharmaceutical Formulation And The Resulting Risks

2

Extractable Study: Applying Different Extraction Procedures And Different Analytical Technologies

3

Toxicological Assessment: Defining Threshold Levels For The Extracted Compounds

4

Leachable Study: Detection, Identification and Quantitation Of Leachables Within The Formulation



Step 1: Evaluating Interactions And Risks

	Likelihood of interaction between packaging component and dosage form		
Degree of concern associated with Route of Administration	High	Medium	Low
Highest	Inhalation aerosols and solution Injections and injectable suspensions	Sterile powders Injection powders Inhalation powders	
High	Ophthalmic solutions and suspensions Transdermal ointments and patches Nasal aerosols and sprays		
Low	Topical solutions and suspensions Topical and lingual aerosols Oral solutions and suspensions	Topical powders Oral powders	Oral tablets Oral hard capsules Oral soft gelatin capsules

Adapted from Guidance for Industry; *Container Closure Systems for Packaging Human Drug and Biologics*, US Department of Health and Human Services, Food and Drug Administration, Rockville, MD, May 1999

- What contributes to the high-risk in pharmaceutical packaging?
 - Prefilled syringe containing an injectable drug suspension
 - Interacts with multiple components in the packaging material (plastic barrel, rubber plunger, metal needle) with direct delivery to the bloodstream

Step 2: Extraction Procedures

Parenteral and Ophthalmic Drug Products (PODP)

Solvents should cover a wide range of polarity

Vigorous conditions

No sample dissolving solvents

No material deformation

Temp Hot extraction techniques

Solvents should mimic drug product formulation

	Thermal	N-Hexane	Isopropanol	Isopropanol/ Water	Aqueous pH 2.5	Aqueous pH 9.5
Headspace	X	---	---	---	---	---
Reflux	---	X	X	PC/PVC only	---	---
Soxhlet	---	X	X	---	---	---
Sealed Vessel	---	---	---	55°C for 3d	121°C for 1hr	121°C for 1hr
Sonication	---	---	---	---	X	X

Autoclave conditions: (121°C for 1hr)

Solvent Polarity/Drug Product Similarity

Detection of additives in LDPE (Example)

Known Additives	Sonication pH 2.5	Sonication pH 9.5	Sealed Vessel IPA/Water
Irganox 1010	---	---	X
BHT	---	---	X
Erucamide	X	X	X

Sonication successfully detects erucamide, but no other anticipated additives

Solvents with different polarity provide better understanding of the material

PQRI: Threshold and Best Practices for Parenteral and Ophthalmic Drug Product (PODP)



Agilent Technologies

Compliance Road Show E/L 2015

Step 2 : Threshold Levels and Actions

<i>Extractable Level in Component</i>	<i>Assignment Category*</i>
> 100 µg/g	Structure Confirmed
20 - 100 µg/g	Confident
< 20 µg/g	Tentative

**Assignment category:*

Structure confirmed: identification categories A, B (or C), and D (or E) (see Table 2) are positive.

Confident: sufficient data to preclude all but the most closely related structures.

Tentative: data is consistent with a class of molecule only.

<i>Identification Category</i>	<i>Typical Identification Data</i>
A	Mass spectrometric fragmentation behavior
B	Confirmation of molecular weight
C	Confirmation of elemental composition
D	Mass spectrum matches automated library or literature spectrum
E	Mass spectrum and chromatographic retention index match authentic specimen

Ref: ITFG/IPAC-RS Collaboration Response to FDA MDI Guidance on 2001



Step 3: Toxicological Assessment & Thresholds (PQRI guidelines for OINDP)

- Based on Toxicological Thresholds of Concern (**TTC**) levels from Kroes et al. (2004)
<http://foodcontactmaterials.com/links/ttc.pdf>
- **Safety Concern Threshold: (SCT)** 0.15 μg per day, which is defined as the threshold below which an individual leachable would have a dose so low as to present **negligible safety** concerns from carcinogenic and non-carcinogenic toxic effects.
- **Qualification Threshold: (QT)** 5 μg per day: Threshold below which a given leachable is not considered for safety qualification (toxicological assessments) unless the leachable presents **structure-activity relationship (SAR)** concerns.
- **Analytical Evaluation Threshold: (AET)** is determined by consideration of the **SCT** and the specific drug product delivery configuration (number of doses in a Drug Product vs single dose)

D. Norwood, L.M. Nagao, C.L.M. Stults; **J. Pharma Sci and Tech.**, (2013) 67(5), 413-429

AET: How much sensitivity is required ?

- 0.05 µg of a genotoxic contaminant was extracted from a drug container
 - Weight opthalmic solution container = 1 g

$$\text{AET} = \frac{0.05 \mu\text{g/container}}{1\text{g material/container}} = 0.05 \mu\text{g/g container material}$$

- Safety Concern Threshold (SCT) for the contaminant = 0.15 µg /d
- Applied Dose = 3 opthalmic solution containers/day
- Content = 3 mL/container

- Analytical requirements to detect leachables within the formulation

$$\text{Leachables: AET} = \frac{0.15 \mu\text{g/day}}{3 \text{ doses/day}} \times 1 \frac{\text{dose}}{\text{container}} = 0.05 \mu\text{g/container}$$

$$\text{AET} = \frac{0.05 \mu\text{g/day}}{3 \text{ ml/container}} = 0.017 \mu\text{g/mL}$$



Step #4 Leachables Study:

❑ Detection of leachables within the formulation

- Different techniques of sample prep are required
 - Extraction procedures to reduce the matrix content
 - Liquid/Liquid extraction
 - SPE
 - Extraction of solids with different solvents
 - Headspace analysis
- Recovery studies are required for quantitative analysis

Threshold levels for leachables in drug products

- **Reported** above **1 ppm** (corresponding to 1 µg /mL sample solution)
- **Identified** tentatively above **10 ppm**
- **Structure confirmation** at **20 ppm**

<http://www.fda.gov/downloads/AboutFDA/CentersOffices/OfficeofMedicalProductsandTobacco/CDER/UCM301045.pdf>

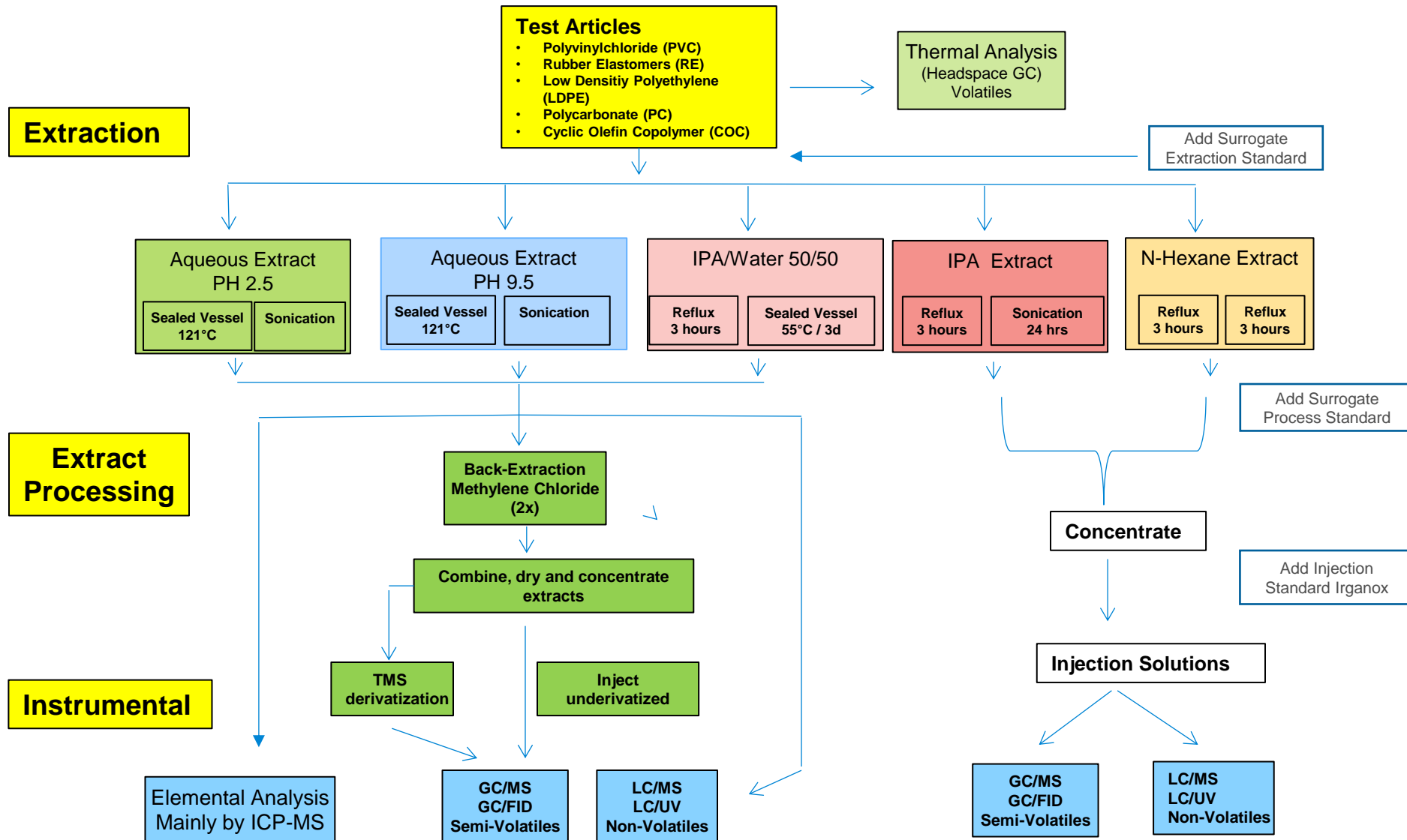


Section #2

- Application Examples: analysis of elemental E&Ls by ICP-MS
- Application Examples: analysis of volatile E&Ls by GC-MS
- Application Examples: analysis of non volatile E&Ls by LC-MS
- Appendix / References



Analytical Workflows: Sample Preparation



Analysis of Inorganic Extractables & Leachables by ICP - MS



Agilent 7800 / 7900 Series ICP-MS



Proposed new ICH and USP methods for elemental impurities: The application of ICP-MS and ICP-OES for pharmaceutical analysis

White paper

Authors

Amir Liba, Ed McCurdy and
Ross Ashdown
Agilent Technologies



Abstract

The United States Pharmacopeial Convention (USP), in parallel with the International Conference on Harmonisation (ICH), is developing new methods for inorganic impurities in pharmaceuticals and their ingredients. The current USP method, <231> "heavy metals limit test", is acknowledged to be inadequate and is due to be replaced with new General Chapters USP <232> (Limits) and <233> (Procedures) in December 2015. The new methods will address the limitations of the current method, extending the list of analytes, reducing maximum permitted exposure limits and taking account of the route of exposure. The new methods will also introduce the use of closed vessel sample digestion and modern instrumental techniques to ensure the accurate recovery and determination of individual analyte concentrations. This White Paper discusses the development of the new USP General Chapters and the ICH Guideline for Elemental Impurities (Q3D) and how Agilent's 7900 ICP-MS and 5100 ICP-OES address the requirements of the proposed new methods.

 **Agilent Technologies**

pub 5990-9382EN, 2014



Agilent Technologies

Published Application Notes



Determination of Chromium in Gelatin Capsules using an Agilent 7700x ICP-MS

Application note

Pharmaceutical

Authors

Miao Jing, Yingping Ni, Yanping Wang and Zhixu Zhang

Agilent Technologies, China



Introduction

Many medications that are administered orally are enclosed within capsules. Frequently both hard-shelled and soft-shelled capsules are made from edible animal protein (gelatin) which is prepared from various animal by-products such as bone and skin. The *Pharmacopoeia of the People's Republic of China* (2010 version) sets a clear standard for the grade of gelatin that can be used for drug capsule production and requires that pharmaceutical companies only purchase capsules from manufacturers that are licensed [1]. There have been recent reports that some companies in eastern China have been making and selling capsules made from cheaper industrial gelatin prepared from discarded leather [2]. Chromium, which is a known carcinogen, and can be toxic if ingested in large quantities, is used in the leather tanning process. Consequently, 20 to 90 times more Cr is typically found in the leather-derived gelatin than in pharmaceutical/edible grade gelatin. As a result, there is a need for a routine, highly sensitive method to



Agilent Technologies

pub 5991-1531EN, 2012



Validating the Agilent 7700x ICP-MS for the determination of elemental impurities in pharmaceutical ingredients according to draft USP general chapters <232> / <233>

Application note

Pharmaceutical

Authors

Samina Hussain

Exova
USA

Amir Liba and Ed McCurdy
Agilent Technologies
USA



Abstract

The United States Pharmacopeia (USP) is developing new General Chapters relating to the determination of elemental impurities in pharmaceutical products and ingredients. USP<232> defines the analyte limits, while USP<233> defines sample preparation options including closed vessel microwave digestion, and recommends the use of modern instrumentation, such as multi-element ICP-MS and ICP-OES techniques. Analytical equipment qualification under USP<233> is based on performance testing, and includes requirements to demonstrate accuracy, repeatability, and the unequivocal identification of analytes. In this paper we present data to illustrate the successful validation of the Agilent 7700x ICP-MS for the measurement of elemental impurities in gelatine capsule samples, according to USP<232> / <233>.

pub 5990-9365EN, 2015



Agilent Technologies

Analysis of Volatile & Semi-Volatile E&Ls by GC-MS



5977B with 7890B GC



7697A Headspace Sampler



7200 Q-TOF



Investigations of Pharmaceutical Products by

- Headspace-GC-MS
- MMI – GC-MS (Multi Mode Injection)

The following pharmaceutical products have been analyzed

- Intravenous (IV) Bag Set & PVC Tubings
- Transdermal Patch
- Liquid Drug Product
- Pressurized Metered-Dose Inhalers (pMDI)



Please read the full story.....

Analysis of Extractable/Leachable Compounds From Plastic Intravenous Bag Sets Using GC/MSD Systems

Application Note

Pharmaceuticals

Abstract

Two Agilent 5977A Series GC/MSD Systems were used for the analysis of extractable and leachable compounds in plastic IV bag sets. Two types of IV bags were investigated: 150-mL dextrose bag (expired) and 1-L sodium chloride bag (warmed). Potentially toxic additives, such as phthalate plasticizers, were shown to have migrated from the IV bag to its infusion solution using the complementation of headspace sampling and liquid injection techniques. High temperature analysis was accomplished using the 7697A Headspace and a 7890A GC coupled with a 5977A MSD. Solvent extracts were analyzed using the MMI 7890A GC coupled with a 5977A MSD. Single ion monitoring (SIM) was used to confirm compound migration.



5991-5616EN

Analysis of Extractable/Leachable Compounds from Transdermal Patches Using GC/MSD Systems

Application Note

Pharmaceutical

Abstract

A lidocaine adhesive patch and film release liner were used to investigate extractable and leachable compounds in transdermal drug delivery systems using two Agilent 5977A Series GC/MSD Systems. Plastic and adhesive additives were identified in acetone, dichloromethane, and hexane extracts using the large volume liquid injection technique. Pharmaceutical ingredients were also identified using high temperature headspace and liquid sampling techniques.

Introduction

Particular interest has been given to extraction techniques in container closure systems (CCS) used in the pharmaceutical industry. Regulators have become increasingly aware of the need to understand whether chemical species can be extracted from the primary packaging material (package with direct contact to the drug product), as well as whether the extracted species (from the package) will appear as leachable species in the drug product. Extractables analysis involves extracting compound from the packaging material using elevated temperatures and solvents related to the packaging composition. Leachables analysis involves identifying compounds in the drug formulation that may have leached from the primary packaging material.

The major source of extractables and leachables are additives that provide physical and protective properties to packaging material, such as flexibility, rigidity, stability, and barrier. Extractables include plastic and elastomeric components, inks and adhesives from coating, and degradation products during processing, storage, and sterilization. Leachables are usually a subset of extractables, however new compounds can form from the interaction between drugs and packaging material.



5991-5605EN

Analysis of Extractable/Leachable Compounds from Generic Liquid Drug Formulations Using GC/MSD Systems

Application Note

Abstract

Pharmaceutical liquid formulations are commonly stored in plastic containers at all risk categories. A pharmaceutical suspension was used as model for investigating compound migration from packaging material. Two Agilent 5977A Series GC/MSD Systems were used. Fatty acid plasticizers were identified using the 7697A Headspace Sampler and a 7890A GC coupled with a 5977A MSD. Phthalate plasticizers were found using the MMI 7890A GC coupled with a 5977A MSD. Single ion monitoring (SIM) confirmed the identification of these plasticizers.



5991-5632EN



Analysis of Extractable Compounds from a Pressurized Metered-Dose Inhaler (pMDI) Using GC/MSD Systems

Application Note

Pharmaceuticals

Abstract

A pressurized metered-dose inhaler (pMDI) is an inhalation device developed for the direct delivery of active pharmaceutical ingredient (API) to the respiratory tract for the treatment of respiratory conditions. Rubber and plastic components in the pMDI are potential sources of extractables by the API/propellant. Therefore, volatile and semivolatile extractable compounds were investigated in these components using two 5977 GC/MSD systems. This application note focuses on identifying extractables in the pMDI using the complementation of headspace GC/MS and MMI GC/MS.

Authors

Diana M. Wong and Roger L. Finor
Agilent Technologies, Inc.
Wilmington, DE, USA

David Weil
Agilent Technologies, Inc.
Schaumburg, IL, USA



5991-6142EN

Agilent Application Note

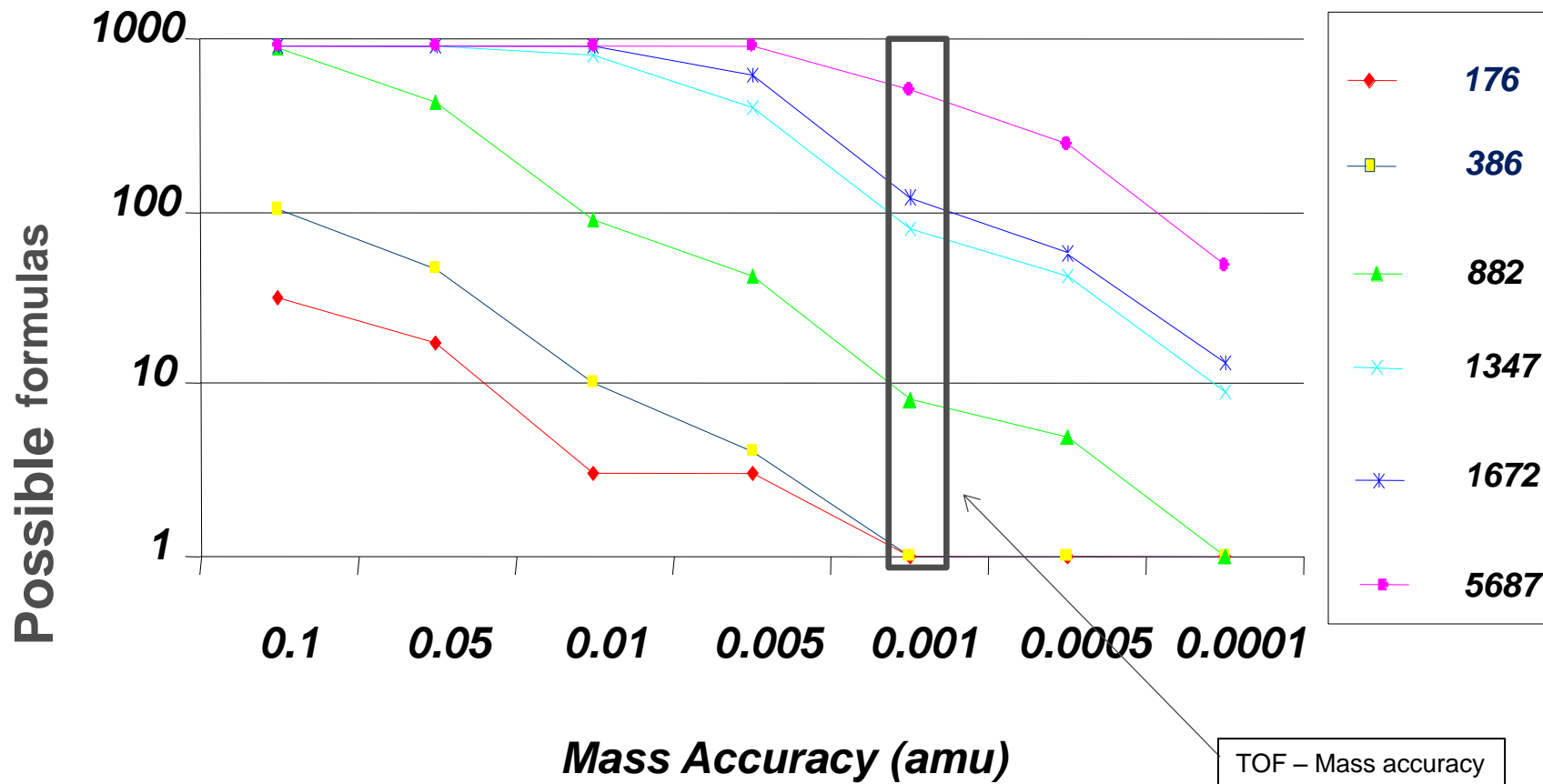
In press.....Agilent Publication Number 5991-6688EN

Differential Analysis in Screening Assays for an E&L Study
Using an Agilent 7200 GC/Q-TOF System Combined with Data
Mining Software

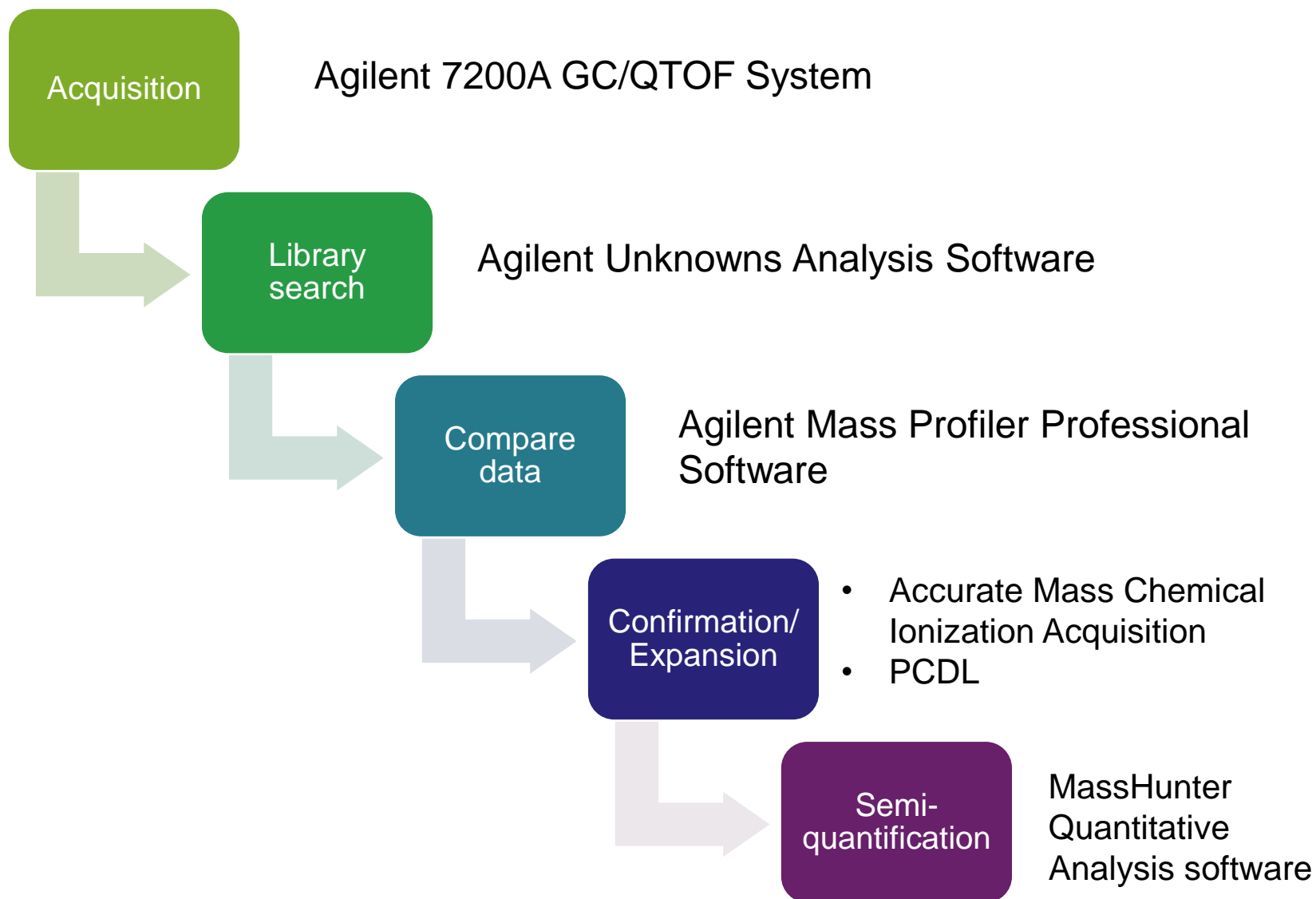


Why TOF Technology?

The Relationship of Mass Accuracy to the Number of Possible Molecular Formulas by Mass

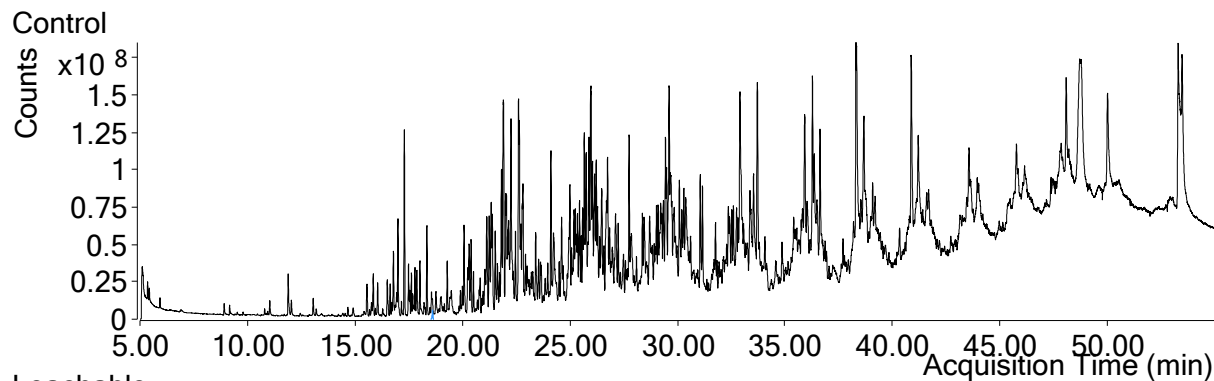


Analytical Workflow

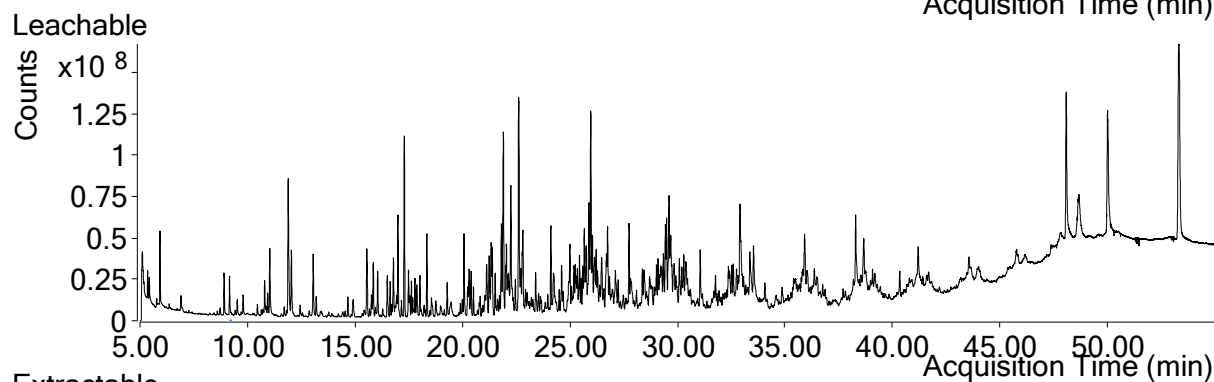


TIC chromatograms in EI mode

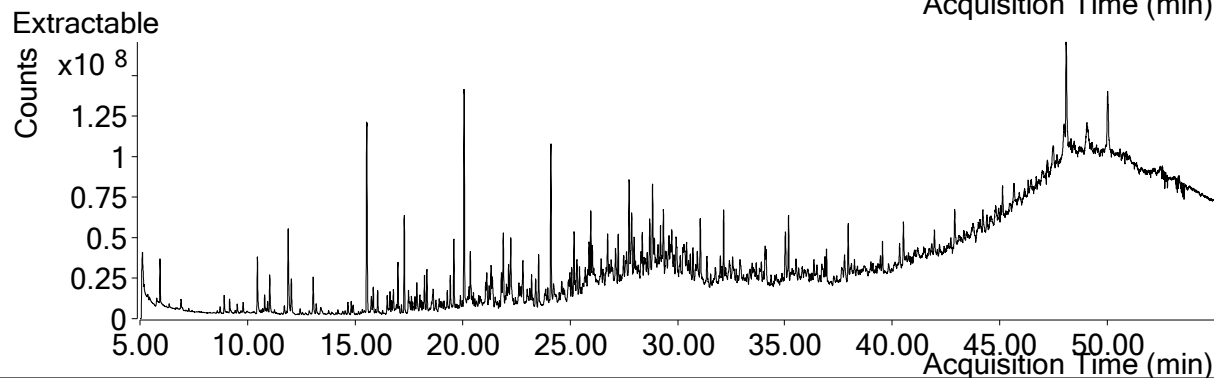
Unheated Formulation
Leachable Sample



Heated Formulation
Leachable Sample

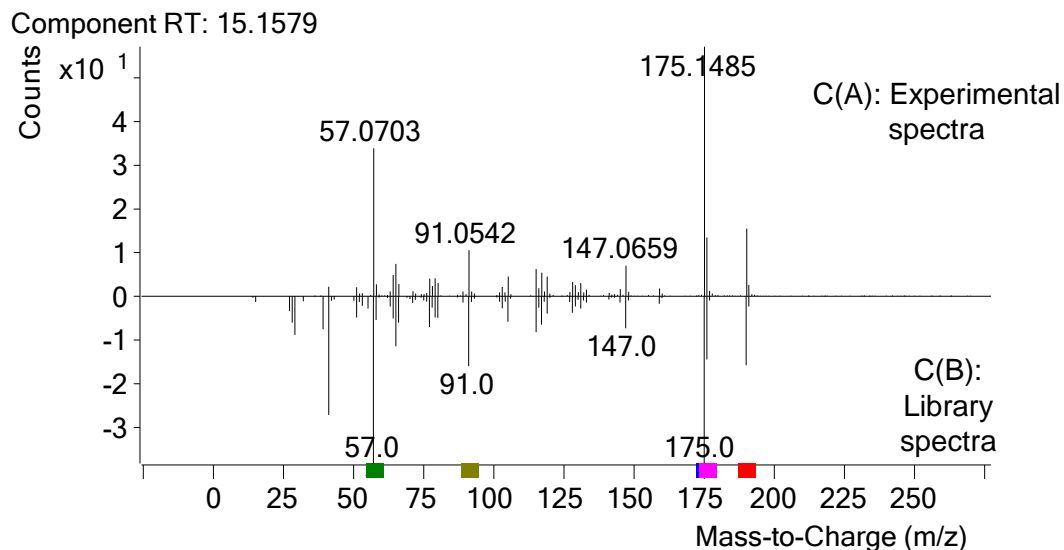
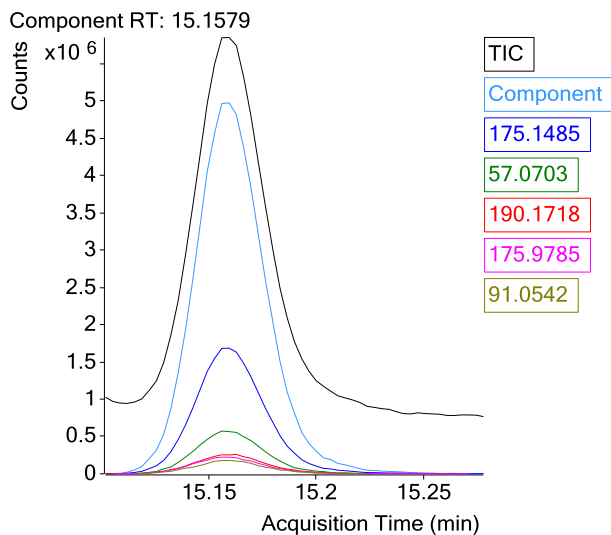
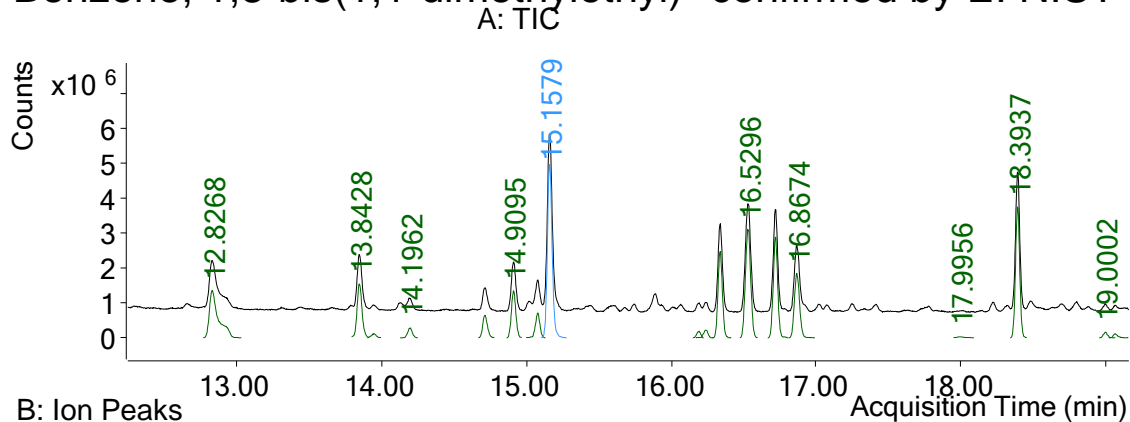


Container Extract
Extractable Sample



Library search by *MassHunter Unknown Analysis Software*

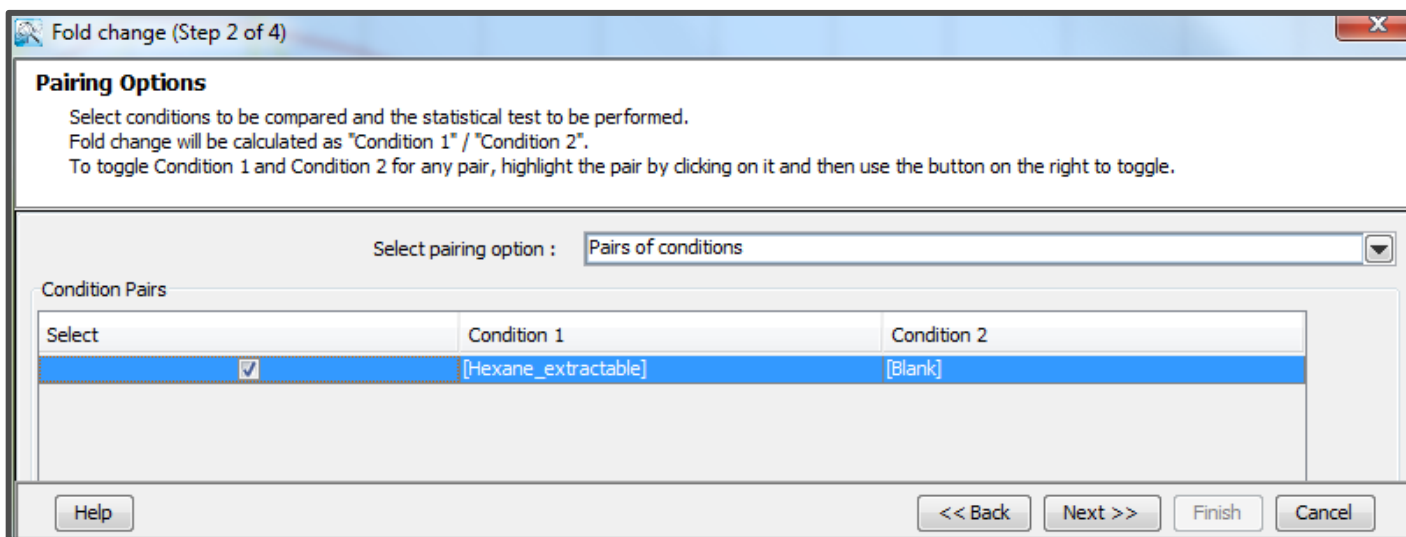
Benzene, 1,3-bis(1,1-dimethylethyl)- confirmed by EI NIST 14.0



Di-Isobutylbenzene is potentially used for polymeric packaging

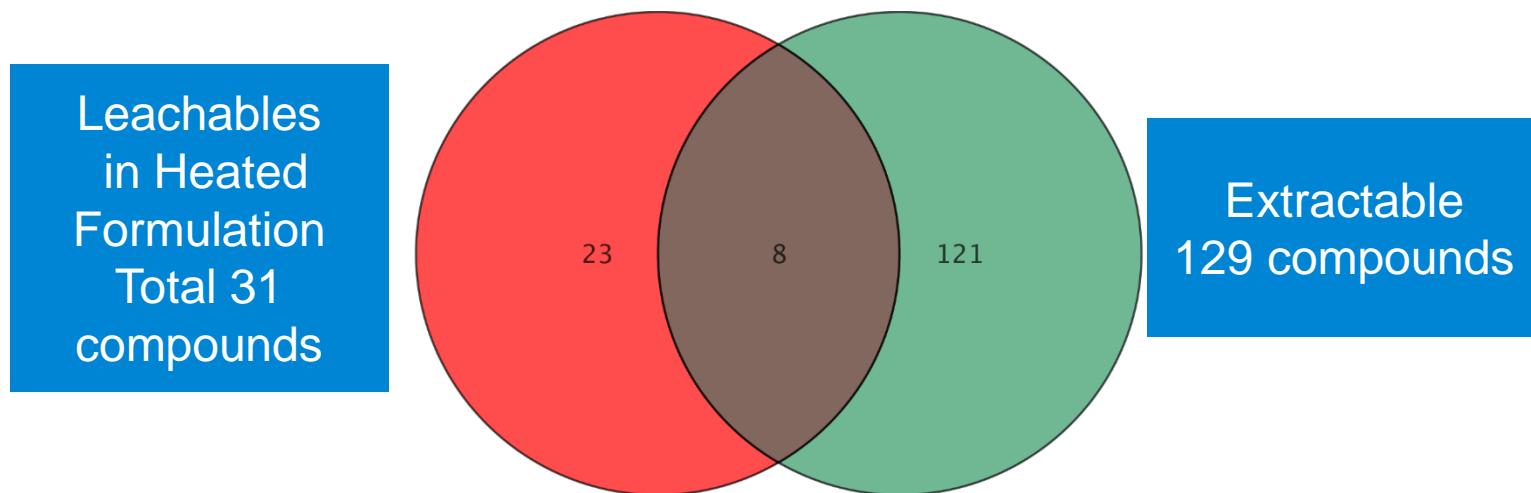
Comparing Datasets to Detect Compounds

Mass Profiler Professional Software



- Traditional blank subtraction could delete an extractable compound by mistake as it might be also present in the solvent, but in lower concentration
- A fold change analysis between the extractable and a blank sample helps to increase the confidence in results

Results: Compounds Found in Different Samples



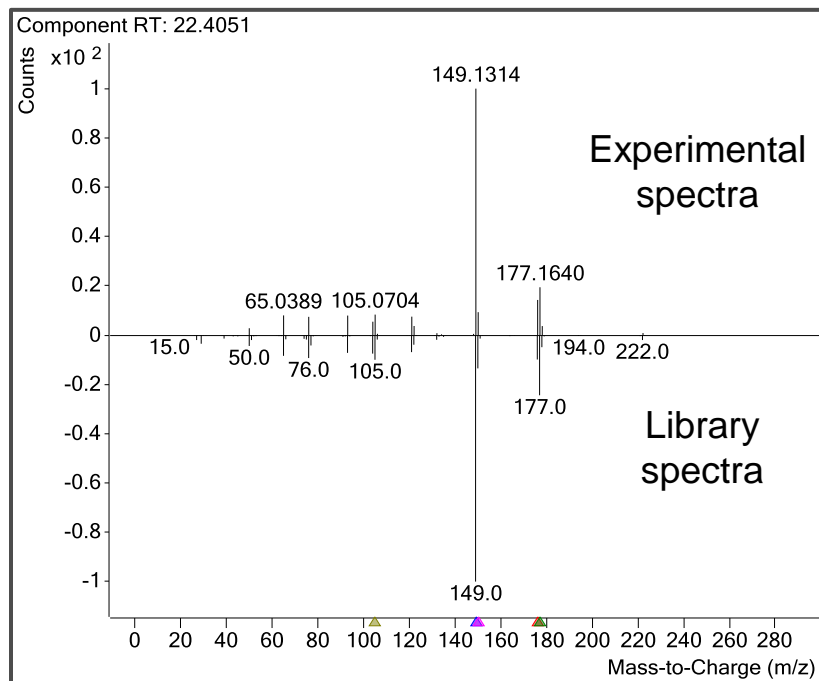
- Benzene, 1,3-bis(1,1-dimethylethyl), and other compounds were also found in the non-heated leachable sample. The origin of these compounds are most likely from the container closure system.

Retention Time	Extractable and leachable compounds	Fold Change in extractable
8.75	Octane, 3,5-dimethyl	UP
15.16	Benzene, 1,3-bis(1,1-dimethylethyl)-	UP
15.75	Dodecane, 4,6-dimethyl	UP
16.19	Tridecane	UP
16.20	Nonadecane	UP
16.87	Cyclohexasiloxane, dodecamethyl-	UP
19.92	Sulfurous acid, pentyl undecyl ester	UP
20.53	Cycloheptasiloxane, tetradecamethyl-	UP



Ambiguity of Results

NIST Library match of diethyl phthalate



Many other compounds also can give the same spectra with matching factor >80

Component RT	Compound Name	Match Factor	Best Hit	Formula	File Name
16.2395	Dodecane, 4,6-dimethyl-	89.8	<input checked="" type="checkbox"/>	C14H30	CE2c.D
15.7427	Dodecane, 4,6-dimethyl-	85.3	<input checked="" type="checkbox"/>	C14H30	CE2c.D
14.7119	Dodecane, 4,6-dimethyl-	86.7	<input checked="" type="checkbox"/>	C14H30	CE2c.D
16.0637	Dodecane, 4,6-dimethyl-	91.0	<input checked="" type="checkbox"/>	C14H30	CE2c.D
13.8455	Dodecane	90.5	<input checked="" type="checkbox"/>	C12H26	CE2c.D
22.4051	Diethyl Phthalate	92.1	<input checked="" type="checkbox"/>	C12H14O4	CE2c.D
28.0820	Dibutyl phthalate			C16H22O4	CE2c.D
28.8504	Dibutyl phthalate			C16H22O4	CE2c.D
8.8540	Cyclotetrasiloxane, octamethyl-			C8H24O4	CE2c.D
12.8274	Cyclopentasiloxane, decamethyl-			C10H30O4	CE2c.D
23.8058	Cyclooctasiloxane, hexadecamethyl-			C16H48O4	CE2c.D
16.8685	Cyclohexasiloxane, dodecamethyl-			C12H36O4	CE2c.D
25.4033	Cyclohexane, undecyl-			C17H34	CE2c.D
23.5054	Cyclohexane, decyl-			C16H32	CE2c.D
20.5230	Cycloheptasiloxane, tetradecamethyl-			C14H42O4	CE2c.D
13.6585	Cyclodecane, methyl-			C11H22	CE2c.D
26.8156	Carbonic acid, octadecyl vinyl ester			C21H40O3	CE2c.D
18.5970	Carbonic acid, eicosyl vinyl ester	86.0	<input checked="" type="checkbox"/>	C23H44O3	CE2c.D

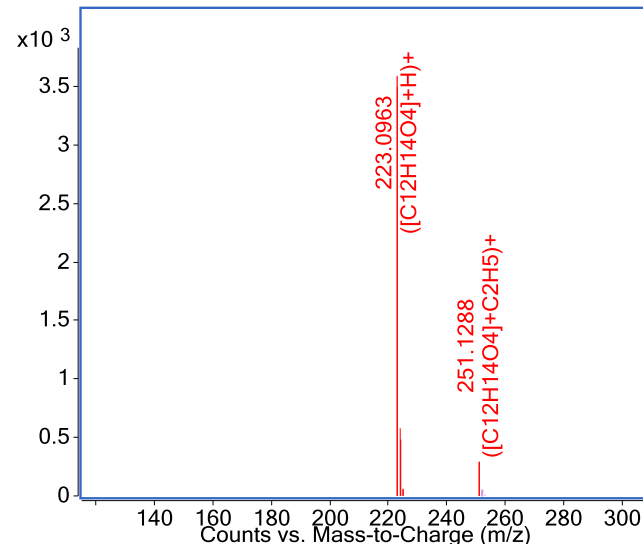
Compound Name	Match Factor	CAS#	Formula
Diethyl Phthalate	93.18	84-66-2	C12H14O4
Diethyl Phthalate	92.71	84-66-2	C12H14O4
Diethyl Phthalate	92.37	84-66-2	C12H14O4
Diethyl Phthalate	91.17	84-66-2	C12H14O4
Diethyl Phthalate	89.81	84-66-2	C12H14O4
Phthalic acid, ethyl isopropyl ester	81.29	1000314-99-6	C13H16O4
Phthalic acid, 2-chloropropyl ethyl ester	80.23	1000356-82-2	C13H15ClO4
Phthalic acid, cyclobutyl ethyl ester	79.43	1000315-41-1	C14H16O4
Phthalic acid, ethyl hex-2-yn-4-yl ester	79.17	1000315-19-0	C16H18O4
Phthalic acid, ethyl pentyl ester	78.97	1000308-93-6	C15H20O4



Eliminating Ambiguity by Chemical Ionization (CI)

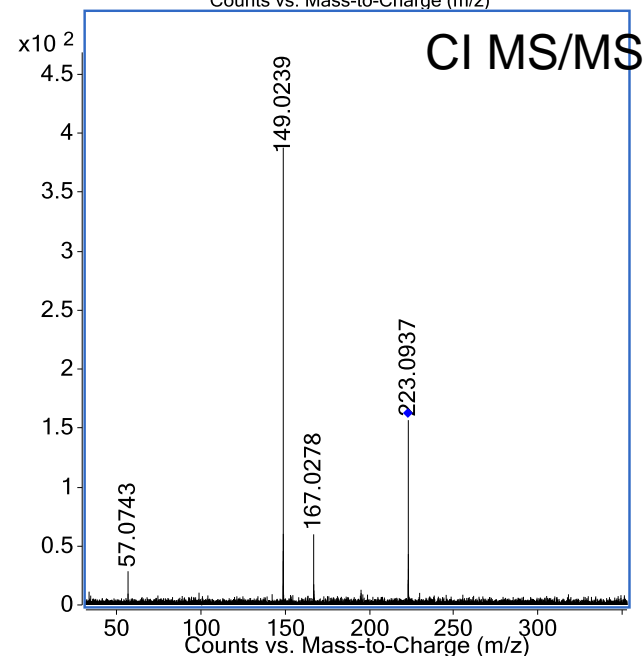
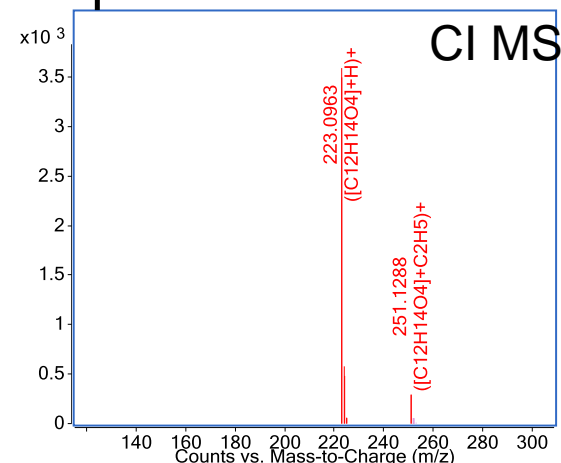
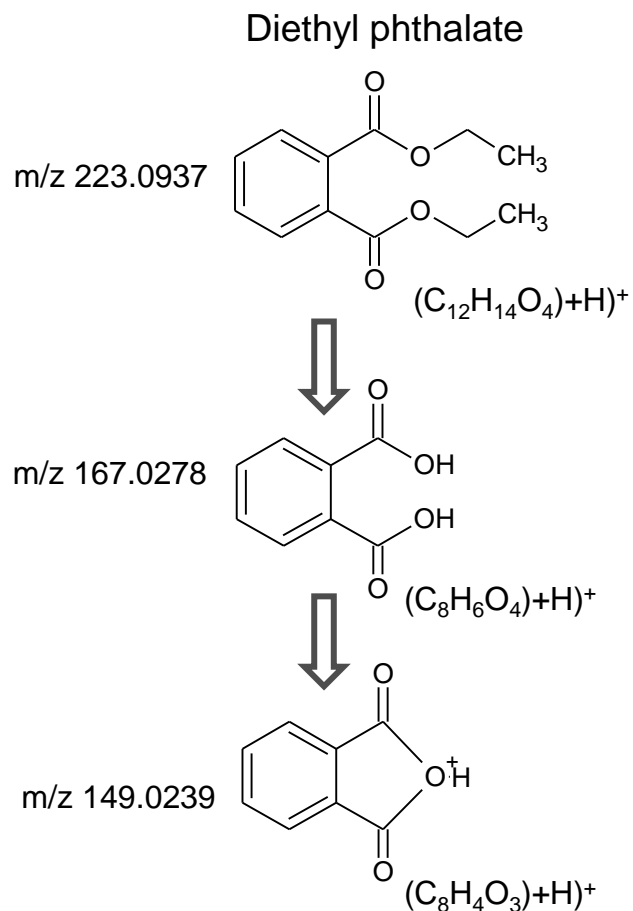
Extractables	Mass	Formula	PPM
.alpha.-Cubebene	204.188	C15 H24	3.52
1,2-Benzenedicarboxylic acid, bis(2-methylpropyl) ester	278.152	C16 H22 O4	4.18
1-Decanol, 2-hexyl-	242.261	C16 H34 O	3.67
2-Methyltetracosane	352.407	C25 H52	-0.6
9H-Fluorene, 9-methylene-	178.078	C14 H10	4.8
Benzene, (1-butylheptyl)-	232.219	C17 H28	3.59
Benzene, (1-butylhexyl)-	218.203	C16 H26	0.07
Benzene, (1-butyloctyl)-	246.235	C18 H30	4.56
Benzene, 1,2,4-trimethyl-	120.094	C9 H12	3.53
Benzene, 1,3-bis(1,1-dimethylethyl)-	190.172	C14 H22	1.44
Benzene, 1,3-dimethyl-	106.078	C8 H10	0.09
Benzene, 1-ethyl-3,5-dimethyl-	134.11	C10 H14	3.55
Benzophenone	182.073	C13 H10 O	3.71
Cyclopentasiloxane, decamethyl-	370.094	C10 H30 O5 Si5	2.85
Diethyl Phthalate	222.089	C12 H14 O4	0.82
Dodecane, 4,6-dimethyl-A	198.235	C14 H30	4.14
Hexadecanal, 2-methyl-	254.261	C17 H34 O	6.11
Naphthalene, 1,6,7-trimethyl-	170.11	C13 H14	5.91
Naphthalene, 2-methyl-	142.078	C11 H10	2.03
(E)-Hex-3-enyl (E)-2-methylbut-2-enoate	182.131	C11 H18 O2	4.99

Diethyl Phthalate (RT 22.4 min)



Eliminating Ambiguity by Collision Induced Dissociation

Fragments & accurate mass for unambiguous compound confirmation



CI MS/MS spectra can be stored in PCDL software to build a custom library.



Creation of libraries from CI-MS/MS data

Database creation

MassHunter PCDL Manager - D:\MassHunter\PCDL\DEP.cdb

File Edit View PCDL Links Help

Find Compounds

Single Search Batch Search Batch Summary Edit Compounds Spectral Search Browse Spectra Edit Spectra

Mass: [M+H]⁺ Neutral [M-H]⁻

Mass tolerance: 10.0 ppm mDa

Retention time: 0.1 min

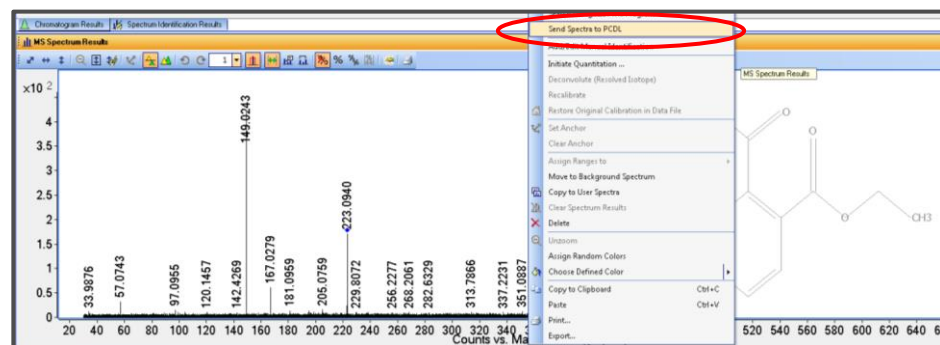
Ion search mode: ☒ Include neutrals ☒ Include anions ☒ Include cations

Formula: Name: Notes: IUPAC: CAS: ChemSpider:

Single Search Results: 1 hit for

Compound Name	Formula	Mass	Anion	Cation	RT (min)	CAS	ChemSpider	IUPAC Name	Spectra
DEP, Diethyl Phthalate	C12H14O4	222.08921	<input type="checkbox"/>	<input type="checkbox"/>	84.662	226	1,2-benzenedicarboxylic acid, 1,2-diethyl ester	0	

Adding the spectra



Library creation

MassHunter PCDL Manager - D:\MassHunter\PCDL\DEP.cdb

File Edit View PCDL Links Help

Find Compounds

Single Search Batch Search Batch Summary Edit Compounds Spectral Search Browse Spectra Edit Spectra

Mass: [M+H]⁺ Neutral [M-H]⁻

Mass tolerance: 10.0 ppm mDa

Retention time: 0.1 min

Ion search mode: ☒ Include neutrals ☒ Include anions ☒ Include cations

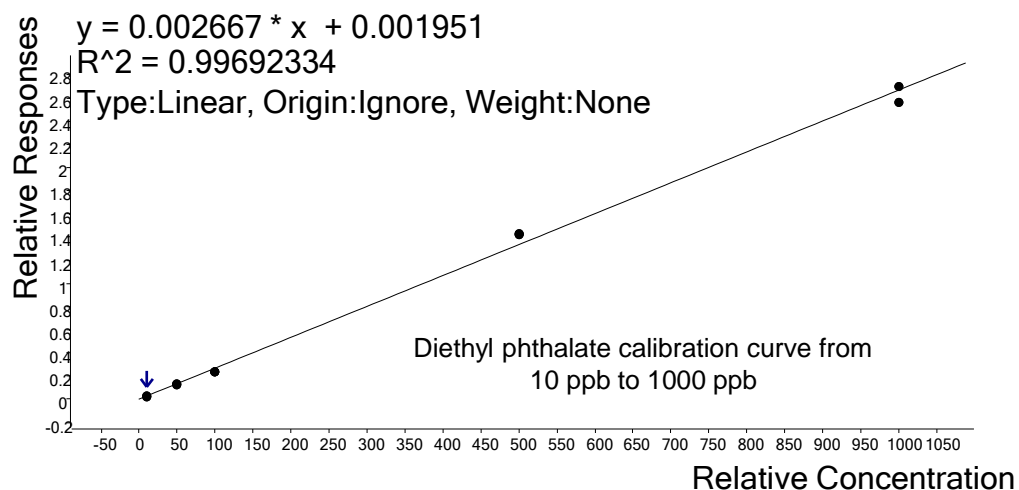
Formula: Name: Notes: IUPAC: CAS: ChemSpider:

Single Search Results: 1 hit for

Compound Name	Formula	Mass	Anion	Cation	RT (min)	CAS	ChemSpider	IUPAC Name	Spectra
DEP, Diethyl Phthalate	C12H14O4	222.08921	<input type="checkbox"/>	<input type="checkbox"/>	84.662	226	1,2-benzenedicarboxylic acid, 1,2-diethyl ester	1	



Semi-quantitative Determination of Impurities



Retention Time	Leachables heated sample	Semi-quantitation estimation (ppb) $\pm 30\%$
6.03	Cyclotrisiloxane, hexamethyl-	88
9.06	Nonane, 2,6-dimethyl-	96
9.12	Sulfurous acid, 2-ethyl hexyl undecyl ester	200
11.42	Octane 5 ethyl 2 methyl	83
13.85	Dodecane	123
15.17	Benzene, 1,3-bis(1,1-dimethylethyl)-	383
16.54	3-Eicosene, (E)-	128
16.79	Tetradecane	58
16.87	Cyclohexasiloxane, dodecamethyl-	127
18.40	Heptadecane, 2,6,10,15-tetramethyl	177
19.31	Dodecane, 2,6,10-trimethyl	44
19.92	Sulfurous acid, pentyl undecyl ester	80
20.75	Phenol, 2,5-bis(1,1-dimethylethyl)-	177
20.98	1-Decanol, 2-hexyl	87
23.89	Heptadecane, 2,6,10,15-tetramethyl	66

- Quantification threshold: 5 ug/day
- Structure confirmed tentatively: 1ppm
- Structure elucidation: 20 ppm
- Semi quantitation: 0.1 ppm to 100 ppm
- The concentration of Benzene, 1,3-bis(1,1-dimethylethyl)- is ~ 0.4 ppm. Based on the daily dosage the consumption (of 9 mL solution/d) is below the quantitation threshold of 5 µg/day.

Analysis of Non-Volatile E&Ls by LC-MS

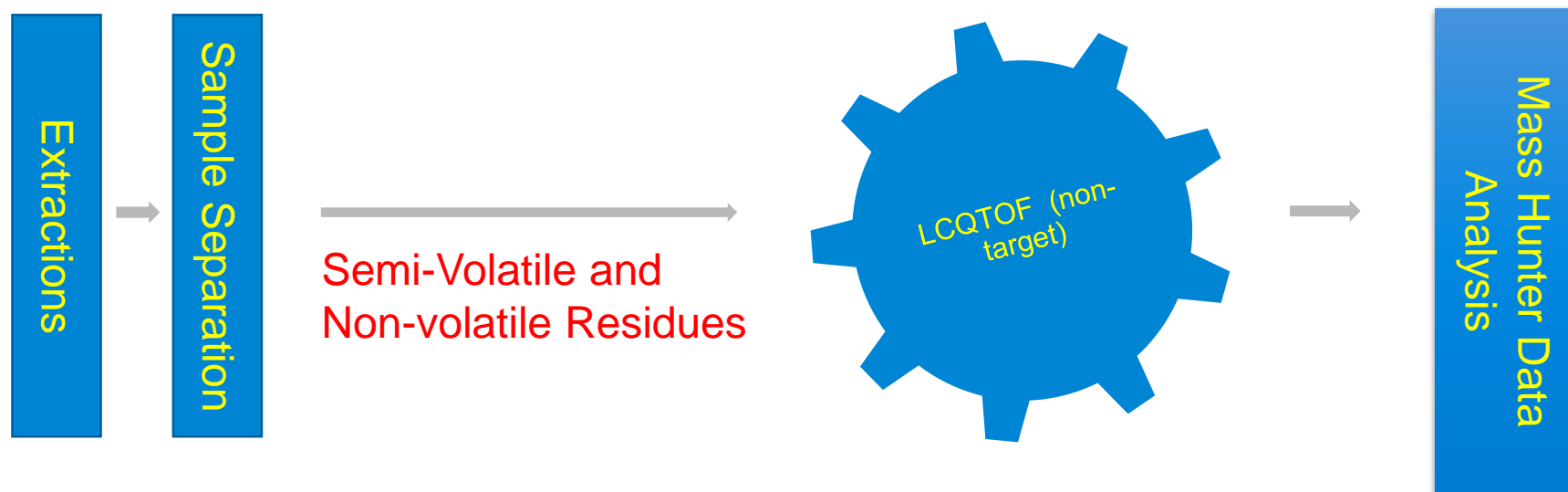
To be published soon.....

Detection and Identification of non volatile E&Ls in an ophthalmic solution by LC-QTOF-MS and MassHunter MassProfiler data mining software

1290 Infinity II UHPLC+ 6500 Series
QTOF System



Extractable/Leachable LC-QTOF Workflow



Samples

- Standards
- Extracts Drug Containers
- Extracts Drug Product

Chromatography

Columns: C18/C8/C3

Organic Mobile Phase:

ACN, MeOH, ACN/IPA,
MeOH/IPA

Varied Buffers: None,
0.1% Formic Acid, 2mM &
4mM NH₄Acetate

Ionization Sources

Jet Stream (ESI),
APCI, Multimode

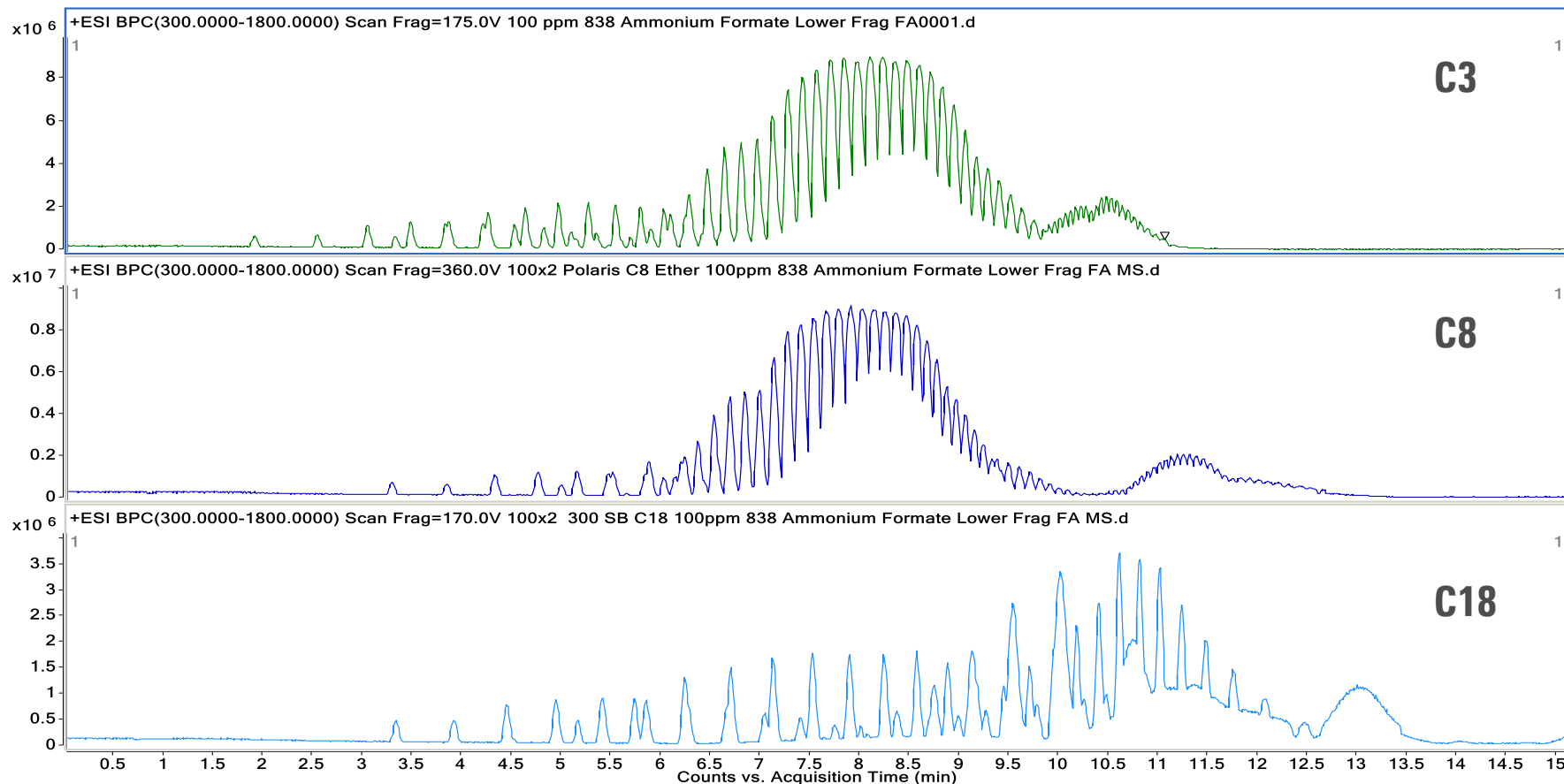
MS-Instrument

6530 / 6545 / 6550



Comparing C3, C8, C18 Separations Using Same Buffers

Base Peak Chromatograms



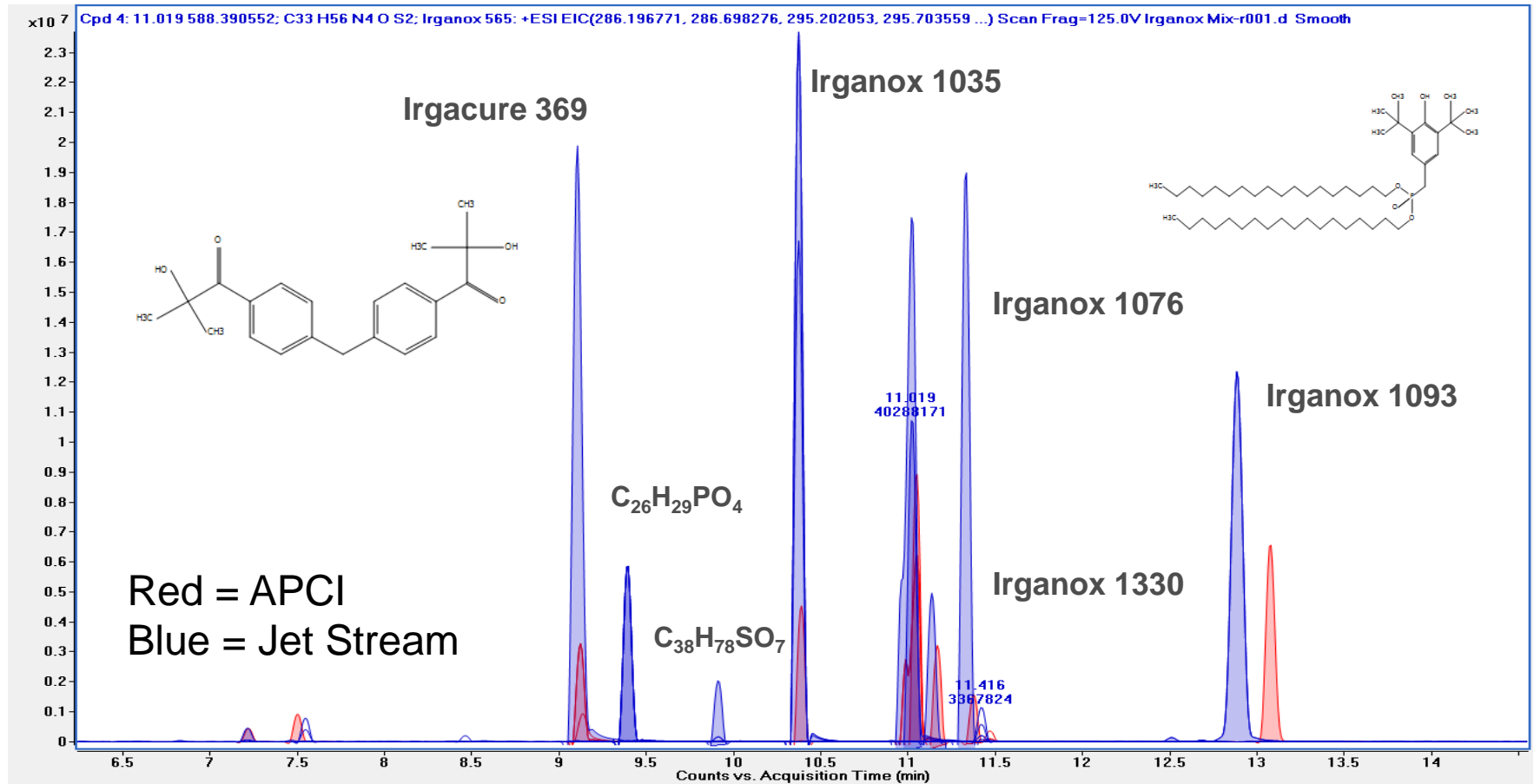
Results: C3 Column Optimum for Higher Mass Extractables
Create orthogonality by using multiple chemistries



Agilent Technologies

Future Directions of Extractable Leachable Analysis David
Weil Oct 2015

Examples: Irganox and Irgacure Mixture



Method developoment for an ophthalmic drug product

System Suitability Test Mix For Method Development

A list of plasticizers has been compiled after literature investigation.
The listed compounds have been selected according to their polarity.

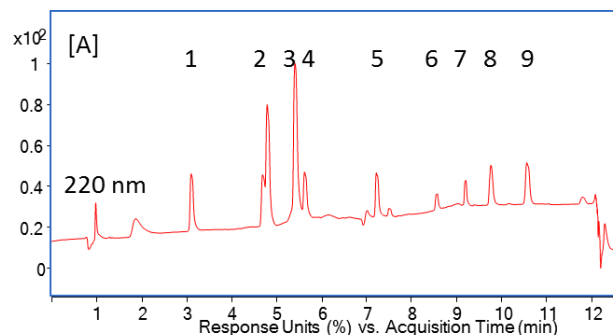
#	Analyte	CAS	Empirical Formula	Monoisotopic Mass
1	Ethyl paraben	120-47-8	$C_9H_{10}O_3$	166.063
2	Irgacure 184	947-19-3	$C_{13}H_{16}O_2$	204.115
3	Irgacure 651	24650-2-8	$C_{16}H_{16}O_3$	256.1099
4	Dipropyl phthalate	131-16-8	$C_{14}H_{18}O_4$	250.1205
5	4-n-Octyl phenol	1806-26-4	$C_{14}H_{22}O$	206.1671
6	Diethyl hexyl phthalate	117-81-7	$C_{24}H_{38}O_4$	390.277
7	Irganox 1010	6683-19-8	$C_{73}H_{108}O_{12}$	1176.7841
8	Irganox 1076	2028-79-3	$C_{35}H_{62}O_3$	530.4699
9	Iragafos 168	31570-04-4	$C_{42}H_{63}O_3P$	646.4515



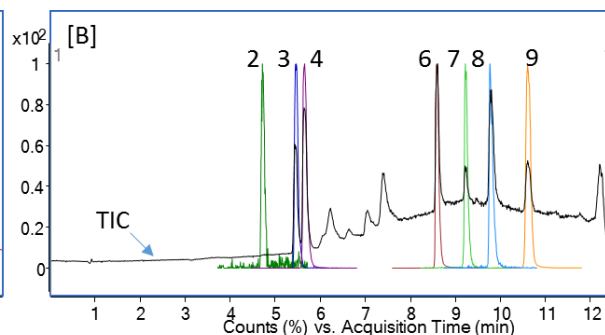
Can I Detect All My Compounds ?

Applying Positive & Negative ES Ionization Mode and UV Absorption

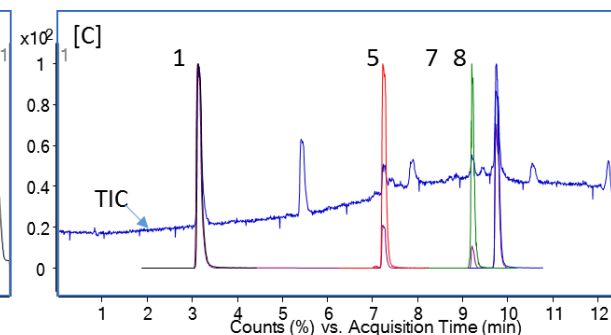
System suitability mix (at 50 ppb level)



UV absorption

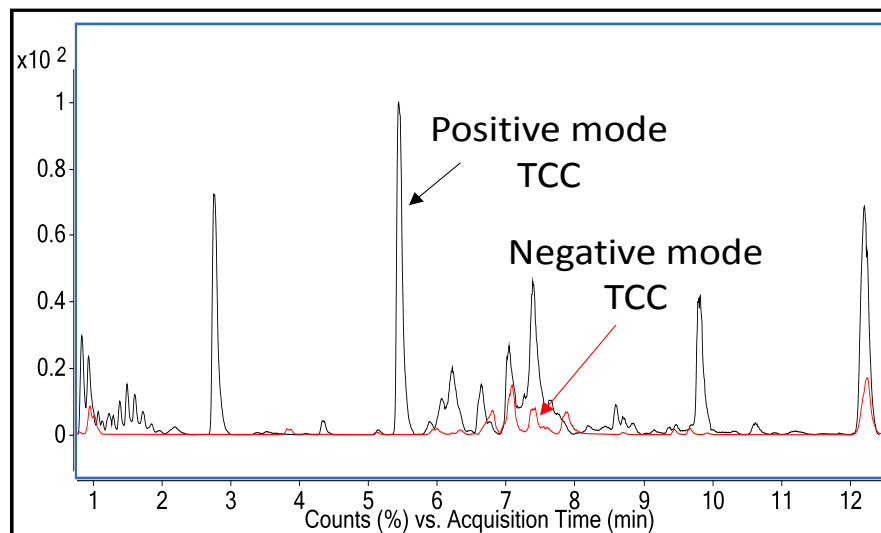


ESI pos; TIC & EIC



ESI neg; TIC & EIC

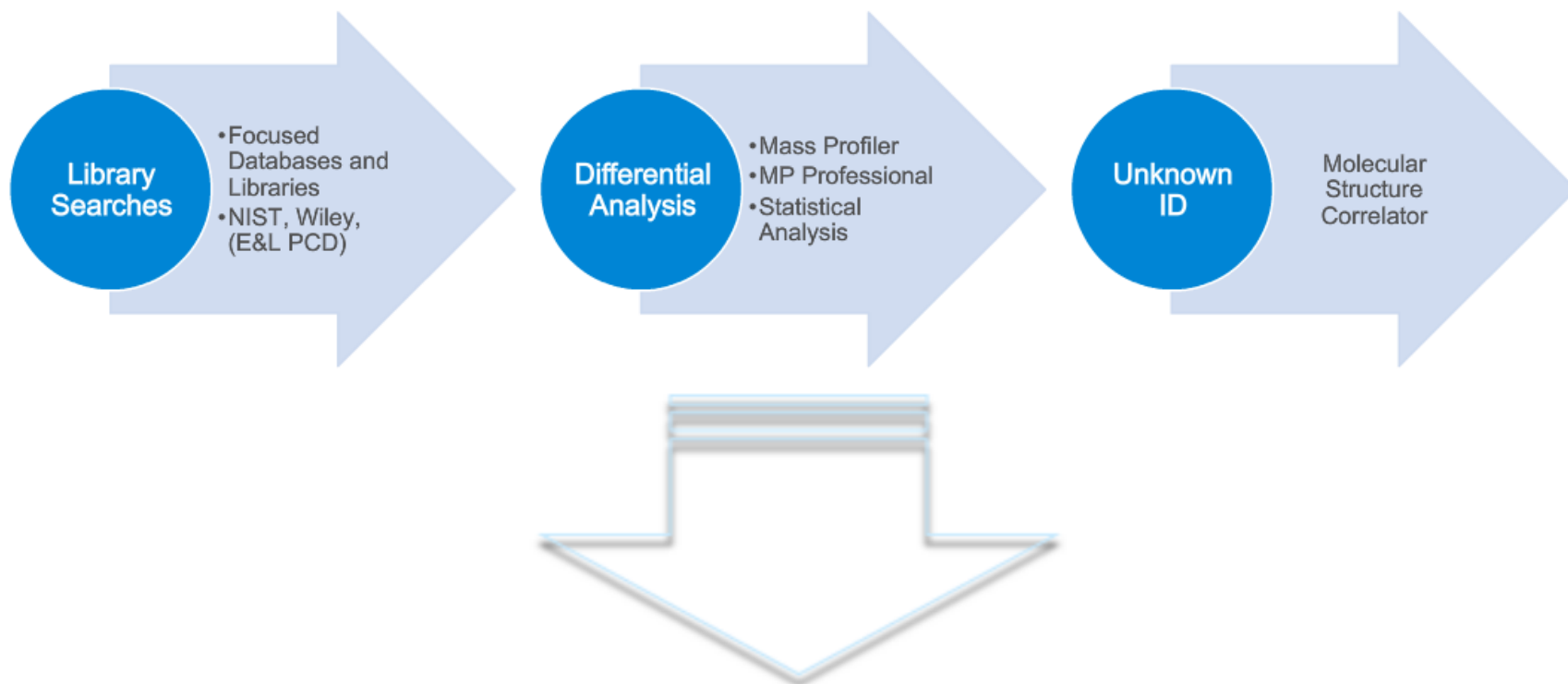
Ophthalmic bottle extract
Overlay of positive & negative TIC
Data acquisition in MS/MS mode



MassHunter Software Tools

Why intelligent data mining software matters....

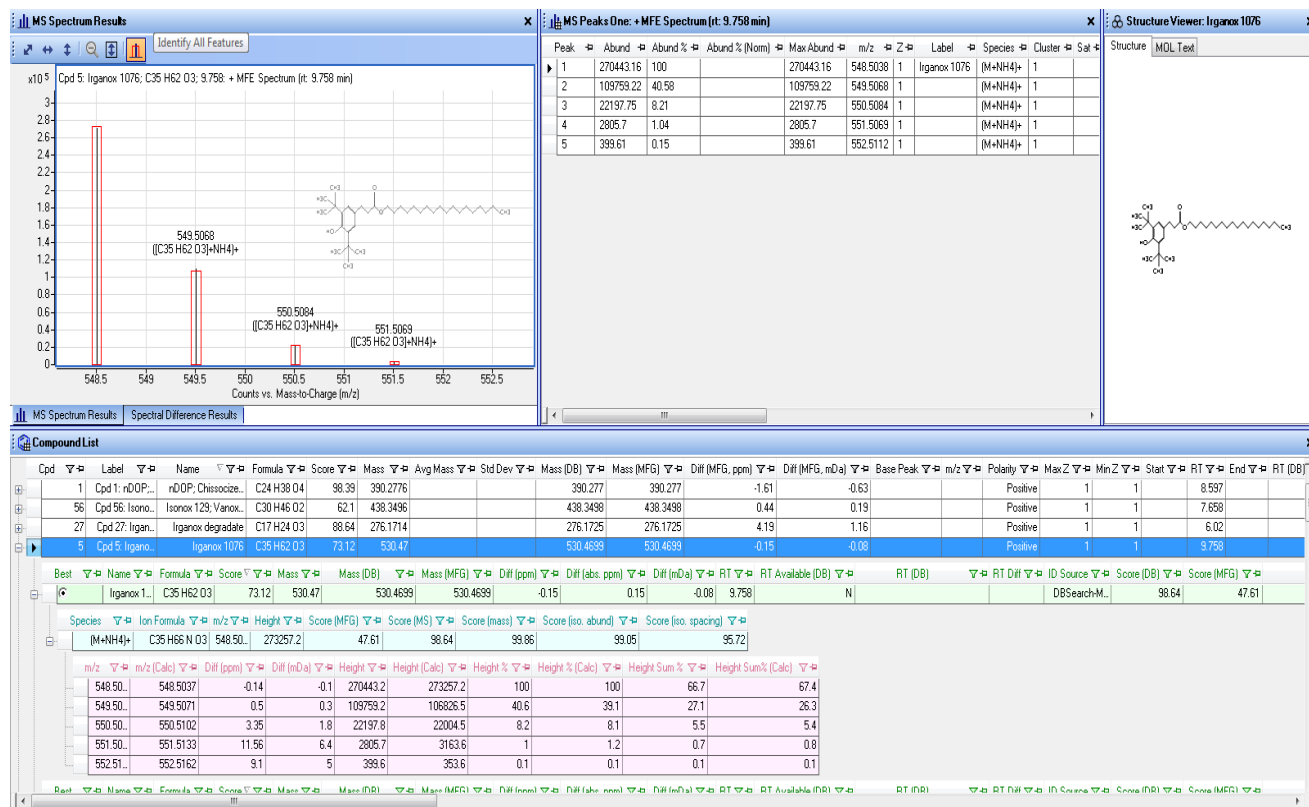
Data mining is an essential step of the analytical workflow and as important as a successful chromatographic separation and detection of organic compounds



MassHunter Quant for Targeted Quantitation



Mass Profiler SW: Database Search & Formula Generation

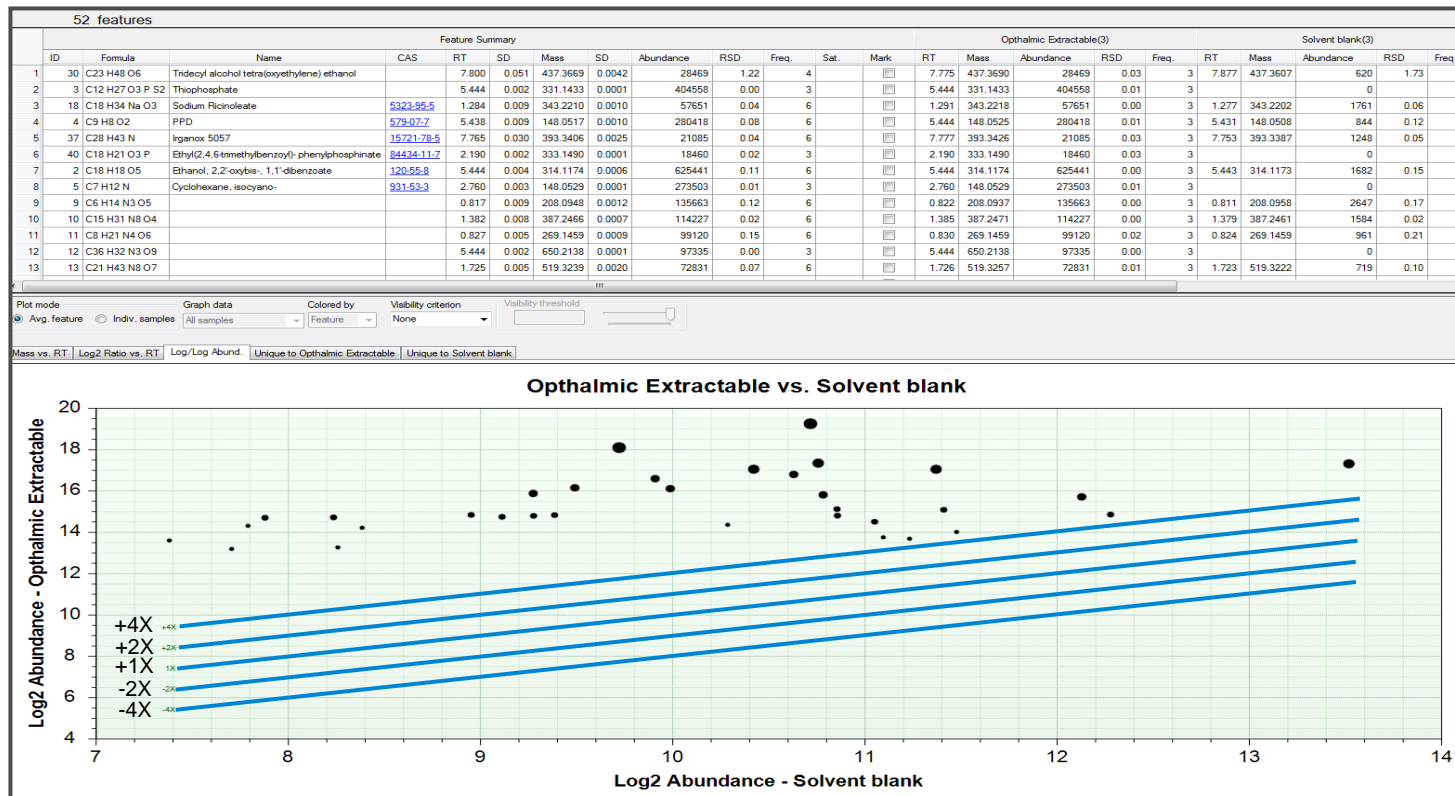


Mass Profiler Software has been used to identify compounds by PCDL comparison
Unknown compounds have been identified by Molecular Structure Correlator Software



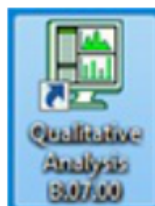
MassHunter Mass Profiler Software (Rev. 7.0)

Differential analysis between the sample and the solvent blank

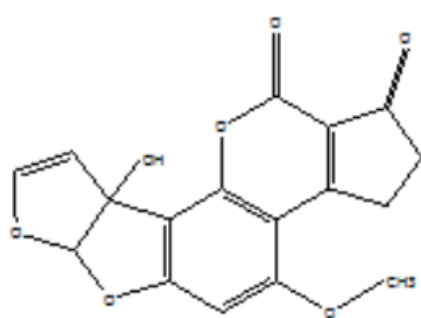


Mass Profiler Software supports the statistical comparison of data sets. After applying a cut off filter of >3,000 and a >4-fold change (abundance by height) and abundance, **66 compounds** (positive and negative ionization modes) have been displayed for the bottle extract

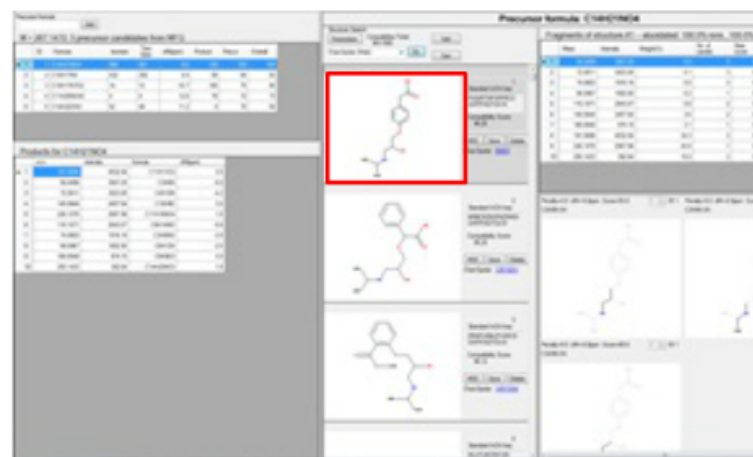
Molecular Structure Correlator (MSC) Workflow



MS/MS spectra in
CEF file



Structures from PCD/L or
ChemSpider, PubChem, etc



Confirmation of structure based
on LC or GC MS/MS data—
**higher confidence
identifications!**

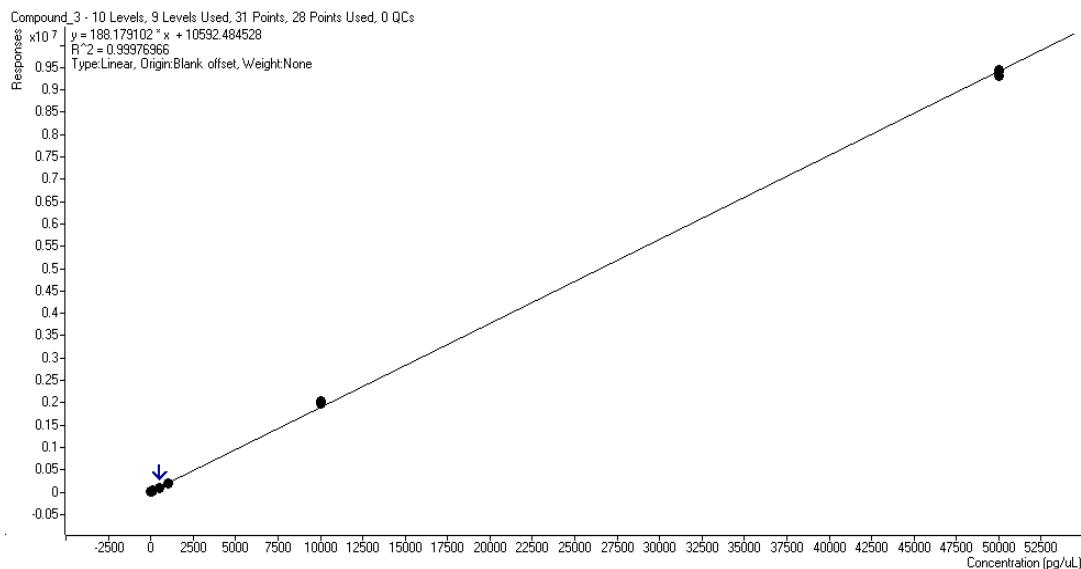


Results: List of identified compounds by PCDL

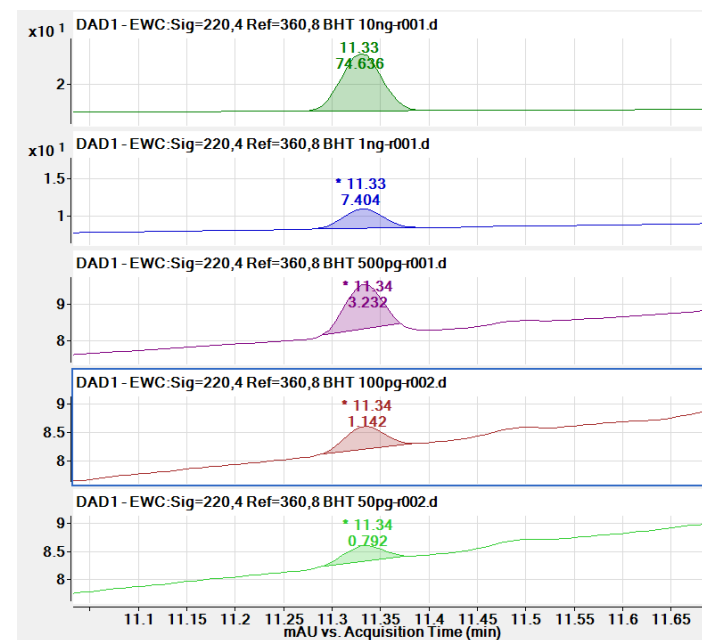
Extractables identified by database	Leachables identified by database
diethylene glycol dibenzoate	Diisononyl phthalate
tridecyl alcohol	Dioctyl phthalate
sodium ricinoleate	Phthalic anhydride
irganox 5057	Methyl-2-benzoylbenzoate
ethyl(2,4,6-trimethylbenzoyl)-phenylphosphinate	Irgacure 907
isocyano cyclohexane	Hexyl Amine
degradant of irganox	Ionox 100
hexadecanoic palmitic acid	Erucamide
Dioctyl Adipate	Glycerol dilaurate
Methyl-2-benzoylbenzoate	Diisodecyl Phthalate
Irgacure 907	Myristyl dimethylamine oxide
Erucamide	Acetic acid, propyl ester
Diisononyl phthalate	
Dioctyl phthalate	
Phthalic anhydride	
Hexyl Amine	
Ionox 100	
Glycerol dilaurate	
Diisodecyl Phthalate	

Semi-quantitative compound determination

Required sensitivity for compound quantification



Quantitation of DEHP $C_{24}H_{38}O_4$
1 pg/μL to 50 ng/μL by Jet Stream ESI



UV Detection of BHT
(Butylated Hydroxy Toluene)
at 220 nm limit around 50 ppb



Results : Semi-Quantification of identified E&Ls

Leachables	ppm $\pm 30\%$
Sansocizer DINP	1.41 ± 0.43
N-DOP	2.48 ± 0.74
Phthalic anhydride	0.14 ± 0.04
Methyl-2-benzoylbenzoate	0.11 ± 0.03
Irgacure 907	0.02 ± 0.005
Hexyl Amine	0.04 ± 0.01
Ionox 100	0.03 ± 0.01
Erucamide	1.68 ± 0.50
Glycerol dilaurate	0.08 ± 0.02
1,2-Benzenedicarboxylic acid, 1,2-bis(8-methylnonyl)ester	0.16 ± 0.05
Myristyl dimethylamine oxide	0.0009 ± 0.0003
Acetic acid, propyl ester	0.10 ± 0.03





Acknowledgements

David Weil	Senior Applications Scientist
Syed Salman Lateef	Pharma Application Scientist
Diana Wong	GC/MS Applications Scientist
Roger Firor	Senior GC/MS Applications Scientist
Anthony Macherone	Senior GC/MS Applications Scientist
Amir Liba	US SPSD AE Manager



Appendix

- Posters published in 2015
- References

Extractable/Leachable Compound Analysis in Pharmaceutical Packaging Using GC/MSD System

ASMS 2015
Poster # ThP 197

67

Posters for ASMS 2015 on E/L

Enhanced Detection of Oligomers, Polymers and Additives Using Ion Mobility Mass Spectrometry and new Data Mining and Differential Analysis Tools

David A. Weil¹; Caroline S. Chu² / ¹Agilent Technologies, Schaumburg, IL; ²Agilent Technologies, Inc., Santa Clara, CA

ASMS 2015
ThP 630

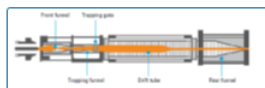


Introduction

The analysis of polymeric/oligomer materials using mass spectrometry has been around since the early 80's using methods such as Field Desorption (FD), Laser Desorption (LD), Matrix-Assisted Laser Desorption Ionization (MALDI) and LC/MS interfaced with a various API methods (APCI, ESI, DESI, ASAP). The complex nature of these materials: (polydisperse, multiple monomers, end groups, degree of branching, MW ranges); challenges determining what stationary and mobile phase, gradients to use. In addition, the polymer mass spectral data with overlapping series of multiply charged species and adducts make data manual data analysis especially challenging. The low electric field ion mobility mass spectrometer (IMS) utilizing separation in the fourth dimension provided a means to simplify the analysis of several classes of structural polymers and also enhanced the detection polymer additives and dyes extracted from a complex mixture. In combination with a new 4D data mining algorithm and differential analysis software to find unique components present in "GOOD" and "BAD" groups of samples. Ion mobility provides a novel means to identify degradation, extractable, leachable impurities in pharmaceutical and consumer products.

Experimental Conditions

The Agilent 6560 uniform low field ion mobility mass spectrometer (drift ion schematic, Figure 2) interfaced with an Agilent 1290 high pressure liquid chromatograph was used for all experiments. Reverse phase separations were carried out using Agilent Stable-Bond-C18 and C18, 2.1 mm x 150 mm, 1.8 µm columns. The mobile phases of water and acetonitrile with 0.1% formic acid was used starting at 1.0% B and ramping to 99% B in 15.0 min with a 5 min final hold time. The flow rates were varied from 0.450 to 1.0 mL/min. The thermally assisted electrospray ionization source operated in positive-ion mode was used for all the experiments. For comparison, data was collected in both QTOF and IM modes.



Structural Information from Drift

Ion mobility mass spectrometry provides insight into the confirmation of metabolites, proteins, glycan, lipids and carbohydrates. For polymers, the adducts used for ionization (H⁺, NH₄⁺, Na⁺, Cs⁺), chain length and degree of branching can lead to confirmation changes as reported by many research groups¹⁻⁴. Star shaped poly(ethylene glycol) PEG polymers are the simplest class of branched polymers with a general structure consisting of several linear chains connected to a central core. These compounds have unique rheological properties that make them useful in drug delivery, thermoplastics and biopolymers. Polymer standards: hydroxyl terminated PEG 2K 4-arm and 10K 8-arm (Figure 1) and Phenyl T and MQ Siloxane resins were analyzed in this study. Confirmation changes were observed as a function of chain length and adducts for these standards.

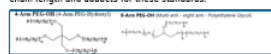


Figure 1: Structures of the 4-arm and 8-arm PEG standards

The extractable/leachable packaging materials extracts samples were prepared using THF solvent. The Beclomethasone dipropionate impurity profiling sample was obtained from a commercial source without additional sample preparation.

Polymer Trend Lines

A major objective in mass spectrometry of oligomeric and polymer materials is determining the chemical compositions of the monomers and end-groups, in addition obtaining information on chain length and molecular weight. Industrial grade materials are rarely "Pure" but typically complex mixtures with monomers, end groups, structures (linear, branched, cyclic). As John McLean's research group⁵ and many others have shown that biomolecules of similar class with follow drift time trend lines. Polymers and oligomers of the same charge state, adduct, end-group composition also follow the trend lines and mass spectral data can be extracted along the trend lines and unique regions using IM Browser software as shown in Figure 3.

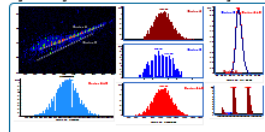


Figure 3: Series A and B from 4-Arm PEG separated by Drift Time and extracted manually along trend line.

Structure/Confirmation Examples

Ion mobility drift heat map 8-arm PEG sample in Figure 4 clearly shows changes in drift time (bends) as a function of molecular weight and chain length. The low molecular weight singly charged linear PEG (shown in top spectrum) clearly separates from the higher molecular weight +3 and +4 charge states branched species (shown on right).

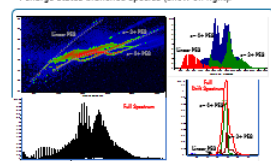


Figure 4: Drift heat map showing change in confirmation as a function of charge state and molecular weight 8-arm PEG sample.

Data Analysis 4D Feature Finding

Manual extraction of mass spectral data (with drift time) is useful but clearly is very time consuming, not automated, may lead to biased reporting and not applicable to applications like proteomics, metabolomics and impurity profiling. A new 4-dimensional (4D) unsupervised data mining algorithm was developed that is based on the 3D Molecular Feature Extraction algorithm, which separates ions as a function of mass/charge, retention time, abundance and drift time⁶. Ions from the same isotope distribution are grouped together and while the adducts are grouped separately with drift time. More details about the algorithm is found in Poster MP168⁷. Results from LC/MS separation of the 4 arm PEG sample is shown in Figure 5.

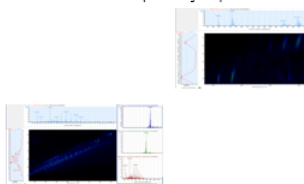


Figure 5: 4D MFE results for LC/MS CE Separation of 4-arm 2K PEG standard

Differential Analysis Mass Profiler

For Extractable Leachable experiments and many others the challenge is to reliably find differences between "Good" and "Bad" lots of sample, at low concentration, with minimum false positive and negative differences. The new 4D molecular feature finding software is integrated into a two sample group comparison with post-alignment filter settings including a feature quality score and an abundance ratio filter.

Either Add Silicone Example showing multiple drift times



Graphical Display Drift Data

Ion Mobility data can be displayed using the Mass Profiler program by: Figure 7 A) m/z versus RT, B) Drift time versus m/z and C) Drift Time versus RT. For 4-Arm PEG we clearly see the presence of minor impurities and drift time changes around RT 6.5 minutes and m/z 600-1000.

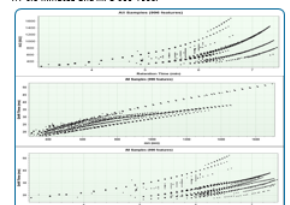


Figure 7A-C: Graphical display 4-arm PEG as a function of Drift time, RT, m/z

Results and Discussion

Extractable Leachable Examples

A major challenge in Extractable Leachable analysis is to identify the presence of low level potentially toxic polymer additives, impurities and dyes in the presence of oligomeric surfactants and polymers. Polymers of similar charge state, adducts, end-groups and composition will follow the same trend lines and mass spectral data can be extracted along those trend lines but additives will be single components and will have different drift times. Mass Profiler was used to identify unique dyes and additives extracted from two different packing materials. A custom accurate mass Extractable Leachable accurate mass database was used to identify copper phthalocyanine C₂₄H₁₈N₄Cu.

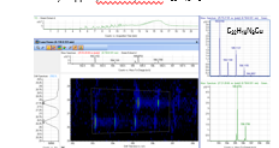


Figure 7: Separation of Possible Copper Phthalocyanine from other extractable using drift and m/z.

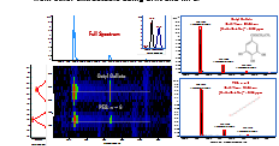


Figure 8: Separation of PEG surfactant and Octyl Gallate antioxidant extractable from expired Claritin medication.

Conclusions

Ion Mobility in combination with high resolution accurate mass, and isotope fidelity using UHPLC separation and data mining with a new 4-D molecular feature extraction (MFE) provides enhanced detection and identification of polymers, oligomers and additives (dyes, antioxidants, pigments) from complex mixtures. Resolved overlapping isotopic patterns.

Impurity Profiling Example

Impurity profiling experiments generally rely on extensive LC method development and/or 2D LC to optimize the separation of low concentration degradation products from the Active Pharmaceutical Ingredient (API). Using Ion Mobility mass spectrometry, two degradation products of Beclomethasone dipropionate (C₂₄H₃₅ClO₆), that overlap chromatographically and have overlapping isotope patterns were separated by drift time. Beclomethasone (C₂₄H₃₅ClO₆) and 21-Dehydro Beclomethasone (C₂₄H₃₃ClO₆) previously not detected using standard LC/MS conditions Figure 9.

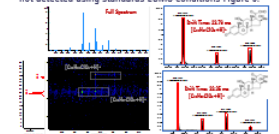


Figure 9: Separation of Beclomethasone (C₂₄H₃₅ClO₆) and 21-Dehydro Beclomethasone (C₂₄H₃₃ClO₆) by drift time

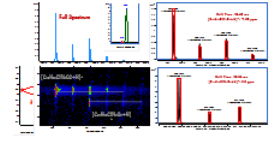


Figure 10: LC/MS Co-elution of degradation products of Claritin C₂₃H₂₇N₂O₂Cl and C₂₃H₂₅N₂O₂Cl

References

- David Clemmer et al. Anal. Chem. 2008, 80, pp 9072-9083
- James A. Solvins et al. Anal. Chem. 2008, 80, 9729-9735
- Scott M. Grayson et al. Macromolecules. 2011, 44, pp 8915-8918
- Edwin De Pauw et al. Anal. Chem. 2014, 86 (19), pp 9892-9700
- John A. McLean et al. Anal. Chem. 2014, 86, pp 2107-2116
- Frank Kuhlmann et al. ASMS 2015 Poster THP 547
- Frank Kuhlmann et al. ASMS 2015 Poster MP 168
- Souvik Kulkarni et al. ASMS 2015 Poster THP 193



Agilent Technologies

Compliance Road Show E/L 2015

Poster for AAPS 2015 on E/L

AAPS 2015
Poster # R6198

A Sensitive Quadrupole Time of Flight Mass Spectrometric Method for Detection and Accurate Identification of Extractables and Leachables

S. Kailasam, S. S. Lateef, S. Khara²

1. Agilent Technologies India Pvt. Ltd, Bangalore, India 2. Agilent Technologies Inc, Santa Clara, USA



Introduction

Drug substances and products can be contaminated by chemicals migrating from primary and secondary packaging materials. Due to the potential impact of these impurities on patient health, the US FDA has issued guidance to the industry on container closure systems for packaging human drugs and biologics. Profiling compounds that can be extracted from the packaging materials, or that have leached (E/L compounds) into a drug substance or product is a complex task due to:

- Wide range of materials used for the construction of primary and secondary containers
- Diversity of physico-chemical properties of the extracted and leached impurities
- Varying concentration of compounds in samples ranging from ng to µg amounts
- Challenges associated with the detection of these compounds in a wide range of different matrices

To overcome some of these analytical challenges, a sensitive high resolution LC/MS method operating in both positive and negative mode has been developed. In addition we have used a statistical approach to detecting and identifying E/L compounds. In this study, a generic method was developed for detecting a set of analytes belonging to different classes of compounds such as plasticizers, photoinitiators, stabilizers, antioxidants, and others. This method was then used for the analysis of E/L compounds in an ophthalmic drug preparation and data processed using a rigorous statistical approach.

Workflow:

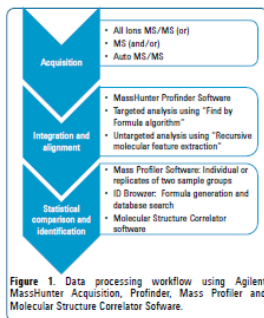


Figure 1. Data processing workflow using Agilent MassHunter Acquisition, Profinder, Mass Profiler and Molecular Structure Correlator Software.

Experimental

Methods

Stock solutions of the standards: ethyl paraben(1), igarcure 18(2), igarcure 65(3), dipropyl phthalate(4), 4,4'-octyl pheno(5), diethyl hexyl phthalate(6), iganox 1010(7), iganox 1076(8) and iganox 18(9), were prepared in isopropanol. A mix of stock solutions diluted in water containing 10 mM ammonium acetate to produce an aqueous mixture containing 1 µg/mL of each analyte. The reverse phase HPLC method and mass spectral parameters were optimized to achieve good chromatographic separation and sensitive analyte detection.

Extractable sample

An ophthalmic medicine bottle, purchased in Bangalore, was washed with water and filled with extraction solvent 1:1 methanol:water. This was incubated in an oven at 55 °C for 72 hours. An aliquot (200 µL) was taken for analysis. The blank used in this study was the extraction solvent.

Spiked sample for method development and system performance test: 10 µL of the standard spike solution (1 µg/mL) was added to 180 µL of the extracted sample.

Leachable/stressed sample

The ophthalmic drug formulation along with its bottle, was heated as is at 60 °C for 24 hrs and analyzed as is. The control in this study was the ophthalmic drug formulation bottle stored at recommended conditions (non-stressed).

Instrumentation

Agilent 1290 Infinity Binary LC System and an Agilent QTOF 6530A System with a Dual Spray Jet Stream Source was used. The LC and MS conditions used in these analyses are shown in Table 1. The source conditions were optimized to enable sensitive detection of extractable and leachable compounds. The LC/MS analysis was performed in both positive and negative modes.

Table 1. LC and MS instrument parameters

LC conditions	
Column	Agilent ZORBAX RRHD Eclipse Plus C18, 3.0 X 100 mm, 1.6 µm (p/n 959768-300)
Column temperature	50 °C
Mobile phase A	100 µg/L Ammonium acetate in water
Mobile phase B	Methanol
Flow rate	0.5 mL/min
Max. Light 60 nm flow cell	220 nm
Gradient	
Time (min)	% Methanol
0	40
8	100
12	100
Stop	11 min
Post	1.5 min
MS conditions	
Ionization mode	Dual Spray A: B: 50
Drying Gas	15 L/min @ 150 °C
Nebulizer pressure	30 Psi
Sheath gas	11 L/min @ 200 °C
Capillary voltage	300 V
Nucleic voltage	300 V
Fragmenter	140 V

Results and Discussion

Analysis of Standards

A spiked sample was used for method development and system suitability test. Figure 2 shows the extracted ion chromatogram from the spiked sample, as detected at UV 220 nm (A), positive ionization mode (B) and negative ionization mode (C). Many standards were found in positive mode, although a few showed better response in negative mode. The LC analysis with UV detection was performed for performance evaluation, however the MS data are discussed due to requirement for analytical specificity and selectivity.

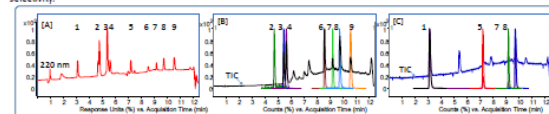


Figure 2. Performance suitability mix analysis using UV at 220 nm (A), accurate mass positive mode (B) and accurate mass negative mode (C). The standards numbering are listed in the method section.

Extractables

The chromatogram from the high resolution LC/MS extractable sample analysis is shown in Figure 3. The total compound chromatogram (TCC) of the sample in both positive and negative modes are also shown. The chemical features in the acquired data were extracted, aligned, and integrated using MassHunter Profinder Software (Figure 4). Any features that failed to integrate can be manually integrated. A total of 175 features from both positive and negative acquisition modes were processed in this manner.

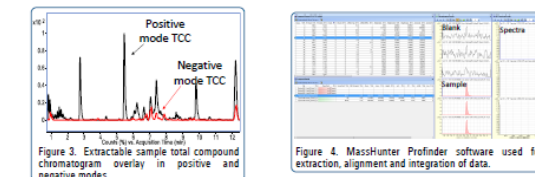


Figure 3. Extractable sample total compound chromatogram overlay in positive and negative modes

Figure 4. MassHunter Profinder software used for extraction, alignment and integration of data.

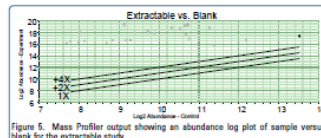


Figure 5. Mass Profiler Software enabled the statistical comparison of data sets. A >4-fold change (abundance by height) and abundance cut off filter of >3,000, displays 66 features (positive and negative ionization modes) for extracted samples

Results and Discussion

Analysis of Leachables

The control and heated ophthalmic formulation samples in leachable study were compared using Mass Profiler software. Statistical comparison using PCA plots is shown in Figure 6. The leachable stressed replicates (red dots) are found to cluster together and distinct from non-stressed replicates (blue dots). These results confirm the significant difference between the non-stressed and stressed leachable samples. Mass Profiler was also used to compare two different sets of samples, the extractables and leachables to determine common and significantly different features (see Figure 7).



Figure 6. Principal Component Analysis of heated (red) and leachable non stressed study samples (blue).

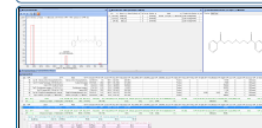


Figure 8. Mass Profiler results showing identified compounds for diethylene glycol dibenzoate.

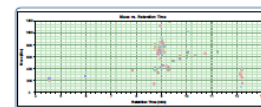


Figure 7. Extractable and leachable sample comparison to identify common compounds and their distribution. A total of 61 compounds were common to both sample types and >2 fold different with each other.



Figure 9. Molecular structure correlator software for confirmation of compounds which were not found in the library.

Identification

The ID Browser functionality within Mass Profiler helped identify compounds using high resolution accurate mass database searching and formula generation from a set of 1,560 compounds (Figure 8). Some of the compounds identified were, triethyl alcohol (oxyethylene) ethanol, sodium ricinoleate, iganox 5057, ethyl(2,6-dimethylbenzoyl)-phenylphosphinate, isocyan cyclohexane, degradant of iganox, and headdecanoic palmitic acid. Library LC/MS-MS spectra matching provided additional compound confirmation. However standards may not be available for many extractable and leachable compounds to build a custom library. Therefore, Agilent Molecular Structure Correlator (MSC) software was used to aid their identification. MSC software proposes possible compound matches by correlating accurate mass/formula of MS/MS fragments with possible in silico fragment ions from proposed structure. Some of the compounds matched with MSC software in leachable sample were: Methyl 2-benzoylbenzoate, Igarcure 907, Eucamide, diethyl adipate, dihydroxyflutranone. MSC helps to provide tentative estimation of the compound identity better than a database search.

Conclusions

- The analysis of a standard mix of nine plasticizers showed that the use of high resolution accurate mass in both positive and negative ionization are required to ensure comprehensive LC/MS analysis of diverse classes of extractable and leachable compounds.
- Integration and alignment verification of samples was performed using Agilent MassHunter Profinder Software, and subsequently samples were compared using Agilent MassHunter Mass Profiler Software.
- The Mass Profiler database search using PCOL and molecular formula generation feature was used to identify compounds. Additional, data analysis of high resolution MS/MS was performed using Molecular Structure Correlator software to confirm the identification of compounds.
- The results of this study show that 66 compounds, including an antioxidant additive, for example, iganox and skin irritants such as: ricinoleate were significantly present at significant concentration in the extractable samples.
- This high resolution and software processing methodology can further be extended to quantification and can also be used for the analysis of samples from drug container closure systems for vendor qualification studies.



Agilent Technologies

Compliance Road Show E/L 2015

Abbreviations

AET =	Analytical Evaluation Threshold
SCT =	Safety Concern Threshold
TDI =	Total Daily Intake
TTC=	Threshold of Toxicological Concern
DP =	Drug Product
OINDP =	Orally Inhaled and Nasal Drug Product
MDI =	Metered Dose Inhaler
QT =	Qualification Threshold
SAR =	Structure-Activity-Relationship

References

- USP Plastic Packaging General Chapters: An Overview, D. Jenke, D. Norwood, Packaging, Storage, and Distribution Expert Committee, USP, http://www.usp.org/sites/default/files/usp_pdf/EN/meetings/workshops/stim_article_661_final.pdf
- USP <1663> ASSESSMENT OF EXTRACTABLES ASSOCIATED WITH PHARMACEUTICAL PACKAGING/DELIVERY SYSTEMS http://www.usp.org/sites/default/files/usp_pdf/EN/meetings/workshops/m7126.pdf
- USP <1664> ASSESSMENT OF DRUG PRODUCT LEACHABLES ASSOCIATED WITH PHARMACEUTICAL PACKAGING/DELIVERY SYSTEMS http://www.usp.org/sites/default/files/usp_pdf/EN/meetings/workshops/m7127.pdf
- Guidelines on Plastic Immediate Packaging Materials, EMEA, European Medicines Agencies Inspections, 2005, http://www.ema.europa.eu/docs/en_GB/document_library/Scientific_guideline/2009/09/WC500003448.pdf
- Extractables and leachables in pharma – A serious issue, F. Moffat, White Paper, http://pat.solvias.com/sites/default/files/whitepaper_extractables_and_leachables.pdf
- Recommendations for Extractables and Leachables Testing, Introduction, Regulatory Issues and Risk Assessment, BioProcess International 5(11):pp36-49 (December 2007), <http://www.bpsalliance.org/wp-content/uploads/2014/06/BPSA-Extractables-101-BPI-Suppl-May-2008.pdf>
- Recommendations for Extractables and Leachables Testing, Executing a Program, BioProcess International 6(1):pp44-53 (January 2008), http://www.bioprocessintl.com/wp-content/uploads/bpi-content/BPI_A_080601AR06_O_76422a.pdf
- The chemical safety assessment process for extractables and leachables associated with packaged pharmaceutical products, D. Jenke, European Pharmaceutical Review, Volume 18, Issue 1, 2013, <http://www.europeanpharmaceuticalreview.com/wp-content/uploads/EPR-Manufacturing-Packaging-Supplement-2013.pdf>
- Metal Leachables in Therapeutic Biologic Products: Origin, Impact and Detection, Shuxia Zhou et al, American Pharmaceutical Review, May 01, 2010, <http://www.americanpharmaceuticalreview.com/Featured-Articles/116570-Metal-Leachables-in-Therapeutic-Biologic-Products-Origin-Impact-and-Detection/>
- Newsletter of the AAPS Aggregation and Biological Relevance Focus Group, May 2011, volume 2, Issue https://www.aaps.org/uploadedFiles/Content/Sections_and_Groups/Focus_Groups/PABCFGnewsMay2011.pdf



References

- HPLC and LC/MS Analysis of Pharmaceutical Container Closure System Leachables and Extractables, D. Norwood et al., Journal of Liquid Chromatography & Related Technologies, 32: 1768-1827, 2009
- Application of the threshold of toxicological concern (TTC) concept to the safety assessment of chemically complex food matrices, M.A.J. Rennen et al., Food and Chemical Toxicology 49.(2011) 933-940
- Leachables and Extractables Handbook, Safety Evaluation, Qualification and Best Practices Applied to Inhalation Drug Products; First Edition, D. Ball, D. Norwood, C. Stults, L. Nagao, John Wiley & Sons, Inc, Published 2012
- Development of Safety Qualification Thresholds and Their Use in Orally Inhaled and Nasal Drug Product Evaluation, Douglas Ball et al., Toxicological Sciences 97 (2), 226 – 236 (2007)
- Regulatory Perspective on Safety Qualification of Extractables and Leachables, Ingrid Markovic, PQRI Workshop, Bethesda (MD), Feb 22, 2011 <http://pqri.org/wp-content/uploads/2015/11/Markovic.pdf>
- Regulatory Perspective on E&L in Biologics: Quality Considerations, Ingrid Markovic, USP/PQRI E&L Workshop, April 28, 2014, Rockville (MD) https://www.usp.org/sites/default/files/usp_pdf/EN/meetings/09_markovich_presentation.pdf
- SAFETY THRESHOLDS AND BEST PRACTICES FOR 6 EXTRACTABLES AND LEACHABLES IN ORALLY INHALED 7 AND NASAL DRUG PRODUCTS, PQRI, 2006, http://pqri.org/wp-content/uploads/2015/08/pdf/LE_Recommendations_to_FDA_09-29-06.pdf
- Current FDA Perspective on Leachable Impurities in Parenteral and Ophthalmic Drug Products, AAPS Workshop on Pharmaceutical Stability, 2011, D. Lewis, <http://www.fda.gov/downloads/AboutFDA/CentersOffices/OfficeofMedicalProductsandTobacco/CDER/UCM301045.pdf>
- Subvisible Particulate Matter, Development in Regulations and Low Volume Methods, Satish K. Singh, AAPS Workshop 2014, http://www.aaps.org/uploadedFiles/Content/Sections_and_Groups/Focus_Groups/Protein_Aggregation_and_Biological_Consequences/PABCFG_Wrkshp20114_Singh.pdf
- Creating a holistic extractable & leachables (E&L) program for biotechnology products, Kim Li, Gary Rogers, Yasser Nashed-Samuel, et al., PDA J Pharm Sci and Tech 2015, 69, 590-619, <http://www.ncbi.nlm.nih.gov/pubmed/26429108>



References

- Perspectives on the PQRI Extractables and Leachables “ safety thresholds and best practices” recommendations for inhalation drug products, D. Norwood, L. Nagao, C. Stults, PDA J Pharm Sci and Tech 2013, 67, 413 – 429
<http://steriletechportal.pda.org/?q=content/pdajpst/67/5/413.full.pdf>
- SAFETY THRESHOLDS AND BEST PRACTICES FOR 6 EXTRACTABLES AND LEACHABLES IN ORALLY INHALED 7 AND NASAL DRUG PRODUCTS, PQRI, 2006, http://pqri.org/wp-content/uploads/2015/08/pdf/LE_Recommendations_to_FDA_09-29-06.pdf
- Current FDA Perspective on Leachable Impurities in Parenteral and Ophthalmic Drug Products, AAPS Workshop on Pharmaceutical Stability, 2011, D. Lewis, <http://www.fda.gov/downloads/AboutFDA/CentersOffices/OfficeofMedicalProductsandTobacco/CDER/UCM301045.pdf>
- Determination of elemental impurities in leachate solutions from syringes using sector field ICP-mass spectrometry, K. Van Hoecke, C. Catry, F. Vanhaecke, Journal of Pharmaceutical and Biomedical Analysis, 77 (2013), 139-144
- Identification and analysis of polymer additives using packed-column supercritical fluid chromatography with APCI mass spectrometric detection, M. Carrot, D. Jones, G. Davidson, Analyst, 1998, 123, 1827-1833
- Analysis of Extractables/Leachable Compounds From Plastic Intravenous Bag Sets Using GC/MSD Systems, D. Wong, R. Firor, Agilent Application Note 5991-5616EN
- Analysis of Extractables/Leachable Compounds from Transdermal Patches Using GC/MSD Systems, D. Wong, R. Firor, Agilent Application Note 5991-5605EN
- Analysis of Extractables/Leachable Compounds From Generic Liquid Drug Formulations Using GC/MSD Systems, D. Wong, R. Firor, Agilent Application Note 5991-5632EN
- Validating the Agilent 7700x/7800 ICP-MS for the determination of elemental impurities in pharmaceutical ingredients according to draft USP general chapters <232>/<233>, S. Hussain, A. Liba, E. McCurdy, Agilent Application Note 5990-9365EN
- Determination of Chromium in Gelatin Capsules using an Agilent 7700x ICP-MS, Agilent Application Note 5991-1531EN
http://www.agilent.com/cs/library/applications/5991-1531EN_AppNote_ICP-MS_7700_pharma_cr_capsules.pdf
- Proposed new ICH and USP methods for elemental impurities: The application of ICP-MS and ICP-OES for pharmaceutical analysis Agilent Application Note 5990-9382EN
- Utilization of Internal Standard Response Factors to Estimate the Concentration of Organic Compounds Leached from Pharmaceutical Packaging Systems and Applications of Such Estimated Concentrations to Safety Assessment, D. Jenke and A. Odufu, Journal of Chromatographic Science, 2012; 50:206-212

