

Comprehensive Extractables Analysis of Medical Grade O-Ring

Dan Ewing,¹ Bill Hurley,² Andrew Feilden,³ Michael Creese,³ Kate Comstock,⁴ and John Schmelzel⁴

¹Parker Hannifin O-Ring Division, Lexington, KY, USA; ²Darcoid Nor-Cal Seal, Oakland, CA;

³Simthers Rapra, Shrewsbury, UK; ⁴Thermo Fisher Scientific, San Jose, CA, USA

Introduction

USP class VI materials are commonly used in medical devices, biopharmaceutical manufacturing process equipment, pharmaceutical packaging, and direct patient-contact applications. For these intended uses, class VI materials need to undergo extractables testing to meet regulatory requirements.



In this study, comprehensive extractables analysis of USP class VI O-rings was conducted using Thermo Scientific™ Q Exactive™ Plus MS and Thermo Scientific™ iCAP™ Q ICP-MS for trace element analysis. Thermo Scientific™ software SIEVE™, Mass Frontier™, and mzCloud™ spectral database were used for data processing.

LCMS Analyses

Sample Preparation

Four different types of medical grade O-rings, A, B, C, and D, were extracted using both water for injection (WFI) and IPA by reflux extraction. For each extraction, 5g of O-ring material and 50 ml of solvent were refluxed for 4 hours. The extracts solutions were analyzed directly by LCMS.

Liquid Chromatography

LC separations were carried out on the Thermo Scientific™ Ultimate™ 3000 RS UHPLC system consisting of: DGP-3000RS pump, WPS-3000RS sampler, TCC-3000RS column compartment and DAD-3000RS UV detector.

Column: Thermo Scientific™ Hypersil™ C18, 2.1x150 mm 1.9 μl Column Temp: 50 °C

LC Mobile phase: A: H₂O/0.1% formic acid B: ACN/0.1% Formic acid

Gradient:

Time (min)	0	3	35	40	40.1	45
B%	10	20	90	90	10	10

Flow rate (μl/min): 400 Injection volume (μl): 10

Mass Spectrometry

MS analyses were carried out on the Thermo Scientific™ Q Exactive™ Plus mass spectrometer using electrospray ionization (ESI) and external calibration. High resolution full scan MS and data dependant top 3 MS/MS data were collected with polarity switching. The resolving power was 70,000 and 17,500 (FWHM m/z200) respectively for full scan and ms/ms. The scan range is 125-1700 amu. HCD stepped normalized collision energy was 25, 35, and 45.



Thermo Scientific™ Q Exactive™ Plus mass spectrometer with a Thermo Scientific™ UltiMate™ 3000 RS UHPLC system

Source Conditions:
 Ionization mode: ESI pos/neg switching
 Sheath gas flow rate: 50
 Auxiliary gas flow rate: 10
 Spray voltage (KV): (+) 3.5 pos / 2.5(-) neg
 Capillary temp (°C): 300
 S-lens RF level: 55.0
 Probe heater temp (°C): 430

Q Exactive Plus MS HRAM Data Acquisition

The High Resolution Accurate Mass (HRAM) data acquisition on Q Exactive Plus MS using full scan/top3 ms/ms with polarity switching generated high quality data, ensuring the detection of structurally diversified compounds at all levels. Figure 1 shows high mass accuracy for pos/neg switching on the fly and the fine isotope pattern of A2. The Δmass1.9954 between A0 and A2 indicate one sulfur present.

FIGURE 1. High Mass Accurate and Fine Isotope Pattern for Pos/Neg Switching

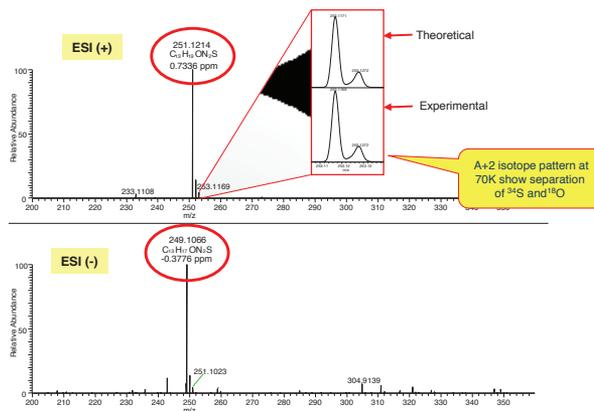


FIGURE 2. Base Peak Chromatographs of Full Scan WFI Extraction (ESI+)

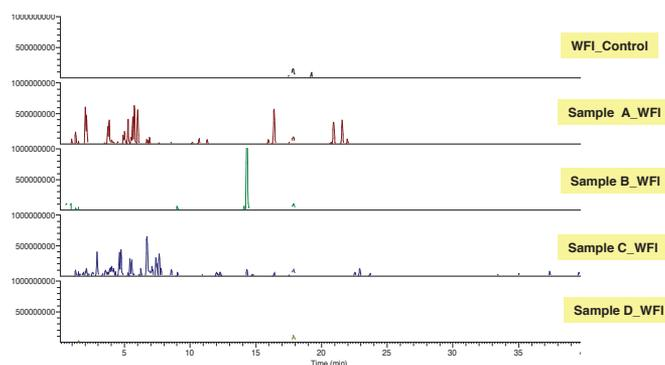
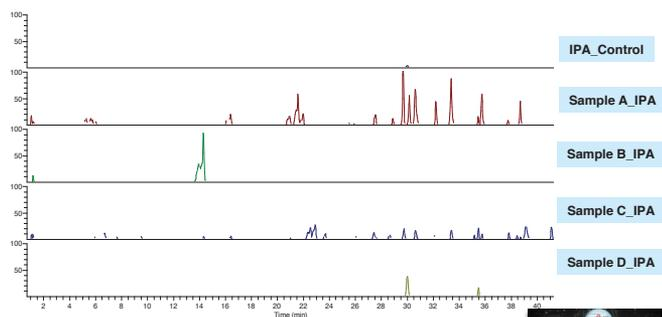


FIGURE 3. Base Peak Chromatographs of Full Scan IPA Extraction (ESI+)



Data Analysis

SIEVE software for Component Extraction and Database Search

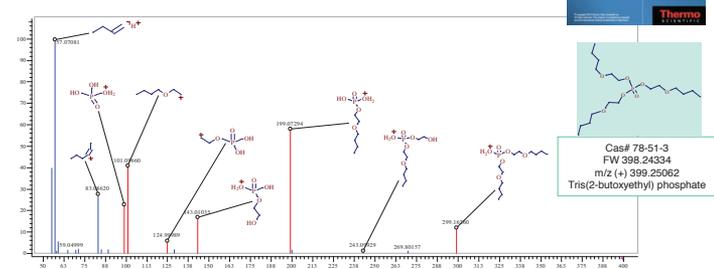
The data was processed using SIEVE 2.2 software for component extraction and differential analysis. The extracted components were filtered using the versatile and customizable filters and then searched against ChemSpider™ database for structure identification.



Component Identification and Structure Elucidation

ChemSpider database searching generated multiple structures for each extracted components. HRAM data with 3 ppm threshold setting reduced the possible structures. To determine the correct structure(s), ms/ms fragments library search was carried out using "Mass Frontier" software, a small molecule structure analysis software. The "HighChem Fragmentation Library" in Mass Frontier software 7.0 has extensive published literature references. For each proposed structure, the "Fragments and Mechanisms" feature in Mass Frontier software was used to generate predicted "fragments and mechanisms" through HighChem Fragmentation Library search, see figure 4. A high degree of correlation between predicted and experimental fragments confirms the proposed structure. Mass Frontier software then automatically annotates the matching fragments based on library search results.

Figure 4. Fragment and Mechanism for Structure Elucidation



mzCloud High Resolution Spectral Database Searching

A search was conducted with "mzCloud". Figure 5 A and B show identification using the ms/ms spectrum search feature.

mzCloud™ is a freely searchable high resolution spectral database at www.mzcloud.org. mzCloud provides several search criteria for small molecule structure identification using tandem mass spectra, including spectra, fragments, precursor ions, etc., all of which can be very useful for unknown structure elucidation.

FIGURE 5-A. Copy MS/MS Spectrum to mzCloud for Substructure Identification

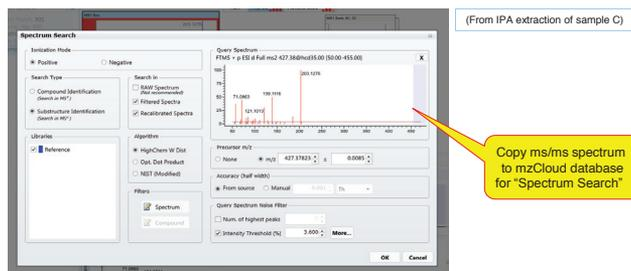


FIGURE 5-B. Spectrum Search Result for Bis (2-ethylhexyl) sebacate

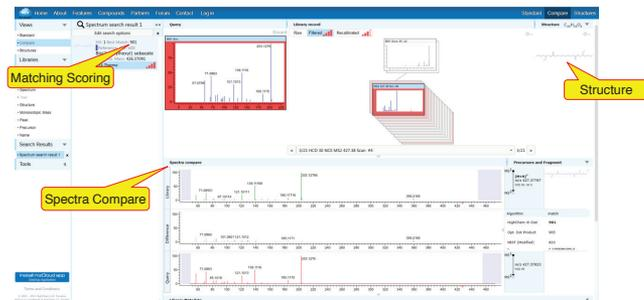


FIGURE 6. Base Peak Chromatogram of WFI Extraction of Sample A

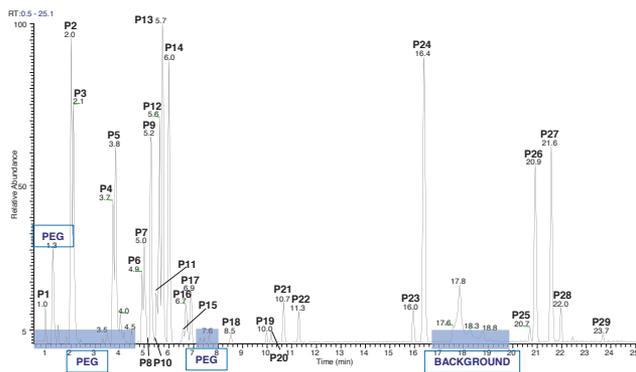


TABLE 1. Components Identified from WFI Reflux of Sample-A (Partial List)

Peak ID	RT	Measured (M+H) ⁺	Calculated (M+H) ⁺	Elemental Composition	Error (PPM)
1	1.0	127.0733	127.0723	C3H6N6	
2	2.0	133.0763	133.0761	C8H8N2	2.1
3	2.1	133.0763	133.0761	C8H8N2	2.1
4	3.7	179.0639	179.0638	C9H10N2S	0.8
5	3.8	179.0639	179.0638	C9H10N2S	0.8
6	4.9	193.0798	193.0794	C10H12N2S	2
7	5.0	193.0798	193.0794	C10H12N2S	1.9
8	5.1	221.0747	221.0743	C11H12ON2S	1.6
9	5.2	237.1058	237.1056	C12H16ON2S	0.7
10	5.3	251.1214	251.1213	C13H18ON2S	0.6
11	5.5	149.0713	149.0709	C8H8ON2	2.1
12	5.6	165.0484	165.0481	C8H8N2S1	1.9
13	5.7	165.0484	165.0481	C8H8N2S1	1.9
14	6.0	251.1214	251.1213	C13H18ON2S	0.7
15	6.6	251.1218	251.1213	C13H18ON2S	1.5
16	6.7	163.1331	163.1329	C8H18O3	0.97
17	6.9	221.1110	221.1107	C12H16N2S	1.3
18	8.5	235.1267	235.1264	C13H18N2S	1.6
19	10.0	265.1372	265.1369	C14H20ON2S	1.1
20	10.2	265.1372	265.1369	C14H20ON2S	1.2
21	10.7	219.1958	219.1955	C12H26O3	1.5
22	11.3	249.1698	249.1697	C12H24O5	1
23	16.0	233.1751	233.1747	C12H25O4	1.5
24	16.4	277.2011	277.2010	C14H28O5	0.2
25	20.7	261.2064	261.2060	C14H28O4	0.4
26	20.9	305.2324	305.2324	C16H32O5	0.1
27	21.6	325.1433	325.1434	C20H20O4	0.1
28	22.0	325.1438	325.1434	C20H20O4	0.3
29	23.7	399.2511	399.2506	C18H40O7P	1.4

TABLE 2. Proposed Structures for WFI Reflex of Sample A (Partial List)

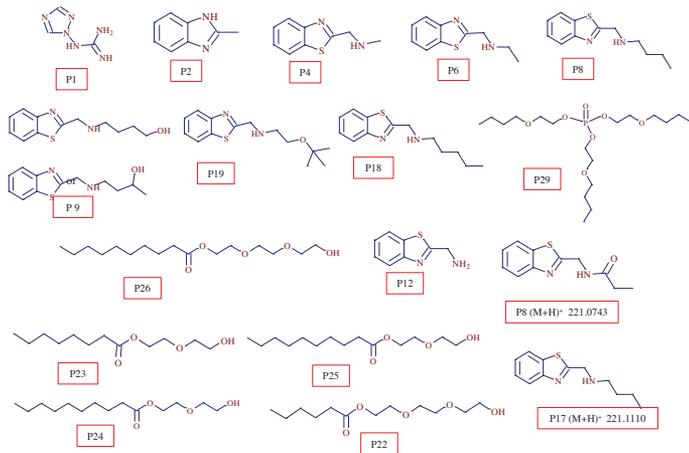


FIGURE 7. ICPMS Results for O-Rings Sample A and B (C & D not shown)

DI Blank	DI EPDM	Nitric Blank	Nitric EPDM	IDL	IDL	IDL	DI Blank	DI Perfluorinated Nitric Blank	Nitric Perfluorinated	IDL	IDL		
6223 (KED)	0.583	241.079	nd	686.431	0.003	0.599	308 (KED)	nd	25.958	0.998	41.121	0.368	1.683
2336a (KED)	1.177	81.969	nd	111.758	0.172	1.724	2336b (KED)	11.377	134.023	nd	38.904	0.132	1.738
4824 (KED)	0.099	48.462	nd	89.891	0.058	0.278	131 (KED)	nd	4.935	nd	35.975	1.359	11.688
396 (KED)	nd	64.572	nd	65.289	0.368	1.051	2340 (KED)	nd	2.462	0.608	14.075	0.060	0.303
2806g (KED)	nd	nd	nd	81.517	0.029	0.288	481 (KED)	nd	1.006	nd	12.762	0.021	0.213
2764 (KED)	nd	nd	nd	68.853	0.000	0.303	569a (KED)	nd	0.334	nd	30.700	0.025	0.147
569a (KED)	nd	nd	nd	38.912	0.015	0.147	569b (KED)	nd	0.334	nd	30.700	0.025	0.147
131P (KED)	nd	3.377	nd	6.442	1.169	11.488	652a (KED)	0.533	11.534	nd	2.728	nd	0.289
630c (KED)	nd	0.789	0.309	2.342	0.004	0.042	652b (KED)	nd	3.346	nd	0.209	0.060	0.599
1378a (KED)	nd	2.684	nd	2.240	0.004	0.035	652c (KED)	nd	3.346	nd	0.209	0.060	0.599
1556a (KED)	nd	0.963	nd	3.893	0.004	0.045	1378b (KED)	nd	45.947	nd	1.599	0.004	0.045
6096 (KED)	nd	0.127	nd	1.902	0.004	0.035	1614a (KED)	nd	1.303	nd	1.000	0.031	0.162
152c (KED)	nd	0.027	nd	1.076	0.003	0.030	1614b (KED)	nd	1.303	nd	1.000	0.031	0.162
6096 (KED)	nd	0.127	nd	1.902	0.004	0.035	909 (KED)	nd	nd	nd	0.240	0.031	0.164
481 (KED)	nd	0.303	nd	0.807	0.021	0.213	909a (KED)	nd	nd	nd	0.240	0.031	0.164
2089b (KED)	nd	nd	nd	0.306	<-0.001	<-0.01	1585a (KED)	nd	0.293	nd	0.234	0.004	0.042
159c (KED)	nd	0.012	0.001	0.200	<-0.001	<-0.01	1585b (KED)	nd	nd	nd	0.234	0.004	0.042
2096 (KED)	nd	nd	nd	0.133	<-0.001	<-0.01	1585c (KED)	nd	0.071	nd	0.189	0.006	0.056
1188a (KED)	nd	nd	nd	0.159	0.006	0.056	1595a (KED)	nd	0.070	nd	0.132	<-0.001	<-0.01
4806a (KED)	nd	nd	nd	0.301	<-0.001	<-0.01	1595b (KED)	nd	0.070	nd	0.132	<-0.001	<-0.01
6101a (KED)	nd	0.049	nd	0.054	0.039	0.199	2089c (KED)	nd	0.046	nd	0.087	<-0.001	<-0.01
1313a (KED)	nd	0.008	nd	0.044	0.004	0.041	1313b (KED)	nd	0.246	nd	0.004	0.041	
9046a (KED)	nd	nd	nd	0.036	0.001	0.014	2098 (KED)	nd	nd	nd	0.028	<-0.001	<-0.01
902c (KED)	nd	nd	nd	0.034	0.002	0.016	79 (KED)	nd	nd	nd	0.074	0.344	0.944
454c (KED)	nd	nd	nd	0.004	0.002	0.042	805 (KED)	nd	nd	nd	0.208	0.004	0.042
74 (KED)	nd	nd	nd	0.076	0.748	118 (KED)	nd	nd	nd	nd	0.423	4.220	2.202
186a (KED)	nd	nd	nd	0.046	0.458	515 (KED)	nd	nd	nd	nd	0.054	0.042	0.245
118 (KED)	nd	nd	nd	0.423	4.226	736a (KED)	nd	nd	nd	nd	0.025	0.225	0.225
515 (KED)	nd	nd	nd	0.025	0.225	736b (KED)	nd	nd	nd	nd	0.025	0.225	0.225
736a (KED)	nd	nd	nd	0.003	0.011	736c (KED)	nd	nd	nd	nd	0.008	0.083	0.083
736b (KED)	nd	nd	nd	0.007	0.072	750a (KED)	nd	nd	nd	nd	0.027	0.072	0.072
750a (KED)	nd	nd	nd	0.027	0.259	750b (KED)	nd	nd	nd	nd	0.027	0.259	0.259
750b (KED)	nd	nd	nd	0.027	0.259	803b (KED)	nd	nd	nd	nd	0.004	0.009	0.009
803a (KED)	nd	nd	nd	0.002	0.019	803c (KED)	nd	nd	nd	nd	0.002	0.019	0.019
1013a (KED)	nd	nd	nd	<-0.001	<-0.01	919a (KED)	nd	nd	nd	nd	0.001	0.007	0.007
1013b (KED)	nd	nd	nd	<-0.001	<-0.01	919b (KED)	nd	nd	nd	nd	0.001	0.007	0.007
1013c (KED)	nd	nd	nd	<-0.001	<-0.01	919c (KED)	nd	nd	nd	nd	0.001	0.007	0.007
1076a (KED)	nd	nd	nd	<-0.001	<-0.01	1076b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1076b (KED)	nd	nd	nd	<-0.001	<-0.01	1076c (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1130a (KED)	nd	nd	nd	<-0.001	<-0.01	1130b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1130b (KED)	nd	nd	nd	<-0.001	<-0.01	1130c (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1130c (KED)	nd	nd	nd	<-0.001	<-0.01	1130d (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1396a (KED)	nd	nd	nd	<-0.001	<-0.01	1396b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1396b (KED)	nd	nd	nd	<-0.001	<-0.01	1396c (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1413a (KED)	nd	nd	nd	<-0.001	<-0.01	1413b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1413b (KED)	nd	nd	nd	<-0.001	<-0.01	1413c (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1475a (KED)	nd	nd	nd	<-0.001	<-0.01	1475b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1475b (KED)	nd	nd	nd	<-0.001	<-0.01	1523a (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1523a (KED)	nd	nd	nd	<-0.001	<-0.01	1523b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1523b (KED)	nd	nd	nd	<-0.001	<-0.01	1575a (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1575a (KED)	nd	nd	nd	<-0.001	<-0.01	1575b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1575b (KED)	nd	nd	nd	<-0.001	<-0.01	1627a (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1627a (KED)	nd	nd	nd	<-0.001	<-0.01	1627b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1627b (KED)	nd	nd	nd	<-0.001	<-0.01	1666 (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1666 (KED)	nd	nd	nd	<-0.001	<-0.01	1679a (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1679a (KED)	nd	nd	nd	<-0.001	<-0.01	1679b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1679b (KED)	nd	nd	nd	<-0.001	<-0.01	1727a (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1727a (KED)	nd	nd	nd	<-0.001	<-0.01	1727b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1727b (KED)	nd	nd	nd	<-0.001	<-0.01	1817a (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1817a (KED)	nd	nd	nd	<-0.001	<-0.01	1817b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1817b (KED)	nd	nd	nd	<-0.001	<-0.01	1817c (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1817c (KED)	nd	nd	nd	<-0.001	<-0.01	1820a (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1820a (KED)	nd	nd	nd	<-0.001	<-0.01	1820b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1820b (KED)	nd	nd	nd	<-0.001	<-0.01	1820c (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1820c (KED)	nd	nd	nd	<-0.001	<-0.01	1939a (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1939a (KED)	nd	nd	nd	<-0.001	<-0.01	1939b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1939b (KED)	nd	nd	nd	<-0.001	<-0.01	1939c (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
1939c (KED)	nd	nd	nd	<-0.001	<-0.01	2027a (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
2027a (KED)	nd	nd	nd	<-0.001	<-0.01	2027b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
2027b (KED)	nd	nd	nd	<-0.001	<-0.01	2027c (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
2027c (KED)	nd	nd	nd	<-0.001	<-0.01	2271a (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
2271a (KED)	nd	nd	nd	<-0.001	<-0.01	2271b (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
2271b (KED)	nd	nd	nd	<-0.001	<-0.01	2288a (KED)	nd	nd	nd	nd	<-0.001	<-0.01	<-0.01
2288a (KED)	nd	nd	nd	<-0.001	<-0.01								



ICPMS Analyses

The samples were prepared by placing the O-rings in 25 ml DI water and 25 ml 2% nitric acid and soaked at RT for 24 hours. The analyses were conducted on Thermo Scientific iCAP Q ICP-MS with He KED (Kinetic Energy Discrimination) interference reduction mode setting. The instrument was standardized at 10 ppb. To determine if trace and potentially toxic metals were leached from the O-rings, the USP <232> Class 1 & 2 elements and additional elements which are commonly analyzed by ICP-MS were determined. The analyses results for the four types of O-rings showed that they are clean of all Class 1 & 2 elements, see Figure 7 for the ICPMS results for sample 1A and B. The system control software Qtegra provides full 21 CFR Part 11 tool set to operate under compliant environments.

Conclusion

This study demonstrated a workflow for rubber O-ring extractable analysis using HRAM data acquisition, data-process software, and database search. The UHPLC/HRAM full MS/HCD MS² with polarity switching data acquisition on Q Exactive Plus MS, coupled with effective process software and database search using SIEVE, Mass Frontier, and mzCloud, significantly increased the confidence and throughput of routine extractable analysis, particularly for unknown components identification and structure characterization.

WFI and IPA extraction profiles of the four types of medical grade O-rings were quickly established by using this workflow. Data not show for sample B, C, and D. The ICPMS results show that the four samples have none or very low level trace element present. GCMS analysis was carried out but the data has not been reported.

References

- ISO10993-5 and -10.
- USP <1663>
- PQRI "L/E Recommendations to the FDA" <http://www.pqri.org/publications/index.asp>

Acknowledgements

The authors would like to thank Parker O-Ring and Darcoide Nor-Cal Seal for providing the O-ring samples, Smithers Rapra for the extractions, and Thermo Fisher Scientific for the MS analyses.

www.thermofisher.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Africa +43 1 333 50 34 0	Denmark +45 70 23 62 60	Japan +81 45 453 9100	Russia/CIS +43 1 333 50 34 0
Australia +61 3 9757 4300	Europe-Other +43 1 333 50 34 0	Korea +82 2 3420 8600	Singapore +65 6289 1190
Austria +43 810 282 206	Finland +358 10 3292 200	Latin America +1 561 688 8700	Spain +34 914 845 965
Belgium +32 53 73 42 41	France +33 1 60 92 48 00	Middle East +43 1 333 50 34 0	Sweden +46 8 556 468 00
Canada +1 800 530 8447	Germany +49 6103 408 1014	Netherlands +31 76 579 55 55	Switzerland +41 61 716 77 00
China 800 810 5118 (free call domestic) 400 650 5118	India +91 22 6742 9494	New Zealand +64 9 980 6700	UK +44 1442 233555
	Italy +39 02 950 591	Norway +46 8 556 468 00	USA +1 800 532 4752

