

# Comprehensive Analysis of Lithium Ion Battery Anode Samples by Ion Chromatography Coupled with High Resolution Mass Spectrometry

Kate Comstock,<sup>1</sup> Rosanne Slingsby,<sup>2</sup> Charanjit Saini,<sup>2</sup> Paul Voelker,<sup>2</sup> Chris Pohl<sup>2</sup>

<sup>1</sup>Thermo Fisher Scientific, San Jose, CA, USA; <sup>2</sup>Thermo Fisher Scientific, Sunnyvale, CA, USA

## Overview

**Purpose:** To demonstrate a workflow using ion chromatography and high resolution mass spectrometry for lithium ion battery (LIB) anode degradation product analysis.

**Results:** LIB anode degradation products were identified from four anode samples..

## Introduction

The lithium ion battery (LIB) is the key component for electric vehicles (EV) and many other electronic devices. The LIB quality directly affects the performance of EV and other devices. Much research has been done in order to improve the performance and increase the efficiency of LIB.

In this study, comprehensive analysis of LIB anode degradation products was conducted using Ion Chromatography (IC) coupled with High Resolution Mass Spectrometry (HRMS).

## Methods

### Sample Preparation

The four LIB anode samples<sup>[1]</sup> were sonicated and rinsed in deionized water. Extracts were filtered through Whatman PP 0.45 µm filters.

### Ion Exchange Chromatography

The ionic separations were carried out on Thermo Scientific™ Dionex™ ICS-2100 IC System using Thermo Scientific™ Dionex™ IonPac™ AG11, AS11 (2 mm) column. Eluent: KOH from 1 to 65 mM in 45 min with gradient. Eluent source: Thermo Scientific™ Dionex™ EGC 500 KOH Cartridge Thermo Scientific™ Dionex™ AERS™ 500 (2 mm) Suppressor<sup>[2]</sup>.

### Mass Spectrometry

The MS analyses were carried out on Thermo Scientific™ Q Exactive™ hybrid quadrupole-Orbitrap mass spectrometer using electrospray ionization in negative mode.

High resolution full-scan MS and top3 data-dependent MS/MS data were collected at resolving power of 70,000 and 35,000 at FWHM m/z 200 respectively. Stepped HCD normalized collision energy (NCE): 30, 45, 60.

## Result

The anode samples were separated by ICS-2100 system based on conductivity, and ions were eluted from the ion-exchange column based on their valences. The eluent was introduced to a Q Exactive mass spectrometer for High Resolution Accurate Mass (HRAM) measurement. The HRAM full scan and ms/ms data provided rich information for confident elemental composition determination and structure characterization. The data was processed using SIEVE for component extraction, followed by ChemSpider and Thermo Scientific HR compound database searching for structure identification. Small molecule structure analysis software (Thermo Scientific™ Mass Frontier™ software) was used to aid with known and unknown structure elucidation.

### High Resolution Accurate Mass Ensured Accurate and Confident Results

Sulfate ( $\text{SO}_4^-$ ) and phosphate ( $\text{H}_2\text{PO}_4^-$ ) have the same unit mass 97.0 amu. High Resolution Accurate Mass (HRAM) data clearly distinguish these two compounds, which ensured unambiguous identification of ion species, especially for unknown degradation product, see Table 1. HRAM MS/MS fragments facilitated structure characterization through accurate fragment ions elemental composition determination, see Figure 2. In addition, HRAM MS/MS readily distinguished co-eluting compounds. See Figure 3.

TABLE 1. Unit mass vs. High Resolution Accurate Mass (HRAM).

| m/z (-)<br>Unit mass | m/z (-)<br>HRAM | Formula (-)                              | Ionic Species   |
|----------------------|-----------------|--|-----------------|
| 97.0                 | 96.9601         | $\text{SO}_4$                            | Sulfate         |
| 97.0                 | 96.9696         | $\text{H}_2\text{PO}_4$                  | Phosphate       |
| 139.0                | 139.0166        | $\text{C}_3\text{H}_5\text{O}_4\text{P}$ | Phosphate Ester |
| 139.0                | 139.0071        | $\text{C}_3\text{H}_7\text{O}_4\text{S}$ | Sulfonate       |

FIGURE 1. Ion Exchange Chromatography and High Resolution Mass Spectrometry Workflow for Lithium Ion Battery Anode Impurity Analysis

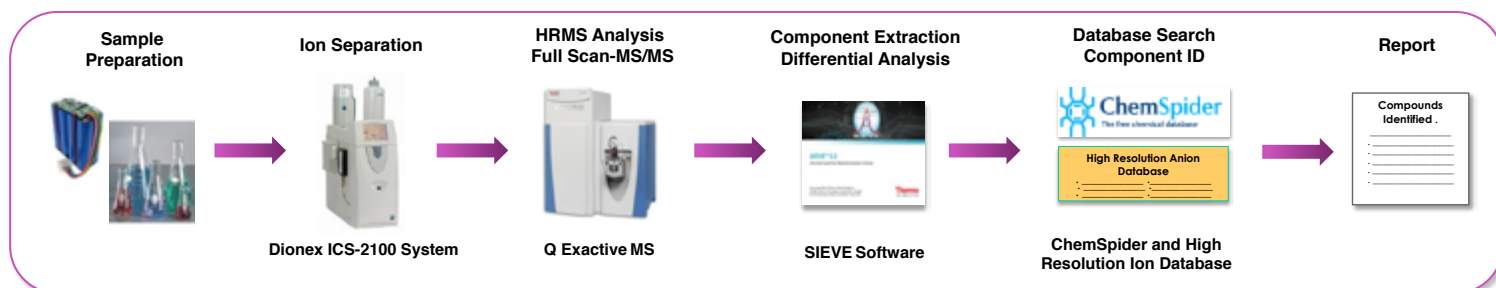


FIGURE 2. HRAM MS/MS Fragments Ensure Confident Structure Characterization

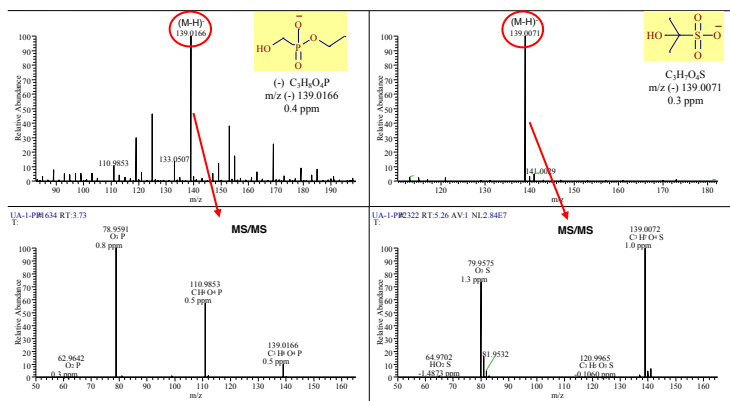
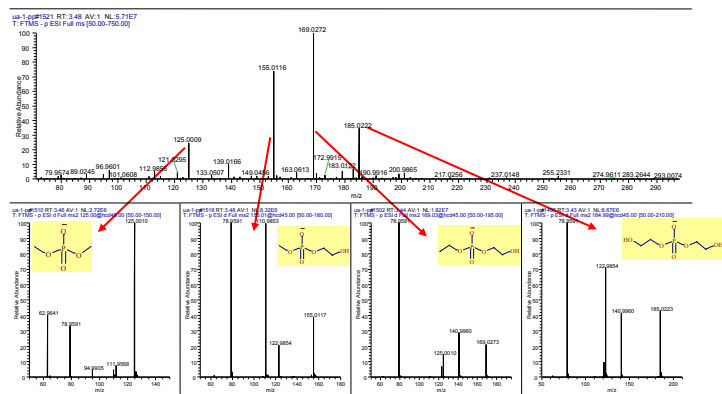


FIGURE 3. HRAM MS/MS Fragments Facilitate Identification of Co-eluting Compounds



HRAM MS/MS Fragments Facilitate Identification of Co-eluting Compounds

The MS negative mode base peak chromatograms of four LIB anode samples with solvent blank and process control are shown in Figure 4 (same scale). Compared with control sample, there were noticeable changes for the other three samples: the peak intensity was either increased or reduced. The data was processed using differential analysis software (Thermo Scientific™ SIEVE™ software) for component extraction and database searching.

FIGURE 4. MS Base Peak Chromatograms (-) of Sample Group

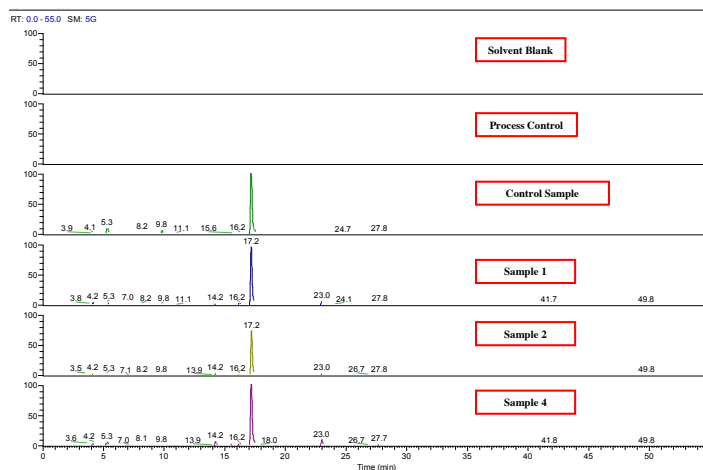
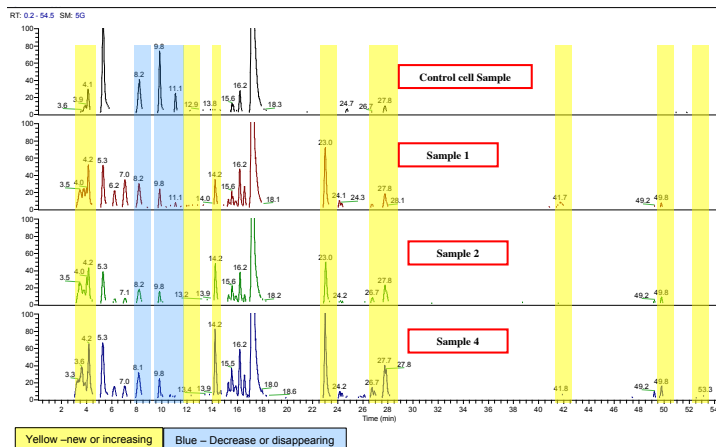


FIGURE 5. Zoomed-in View for Details



SIEVE software results show the extracted components, also the trend intensity and it's details. Figure 6 shows trend intensity of  $m/z$  124.9912 at RT 11.0 min with Elemental Formula  $\text{C}_2\text{H}_5\text{O}_4\text{S}$ .

FIGURE 6. Trend Intensities View for  $m/z$  124.9912 at RT 11.0 minute

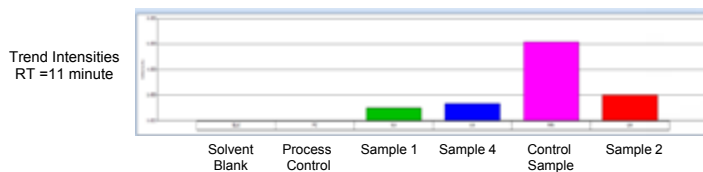
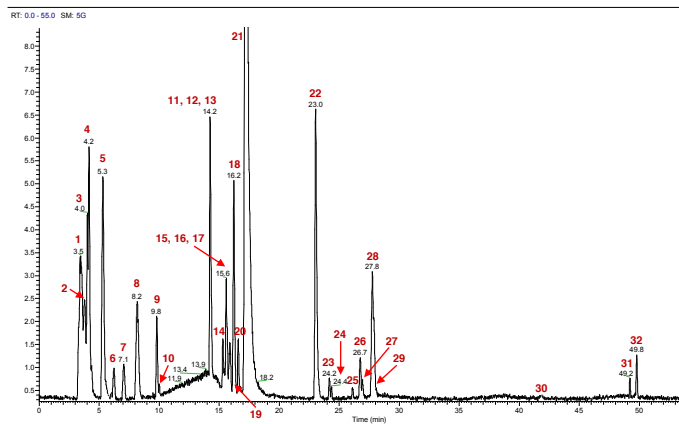


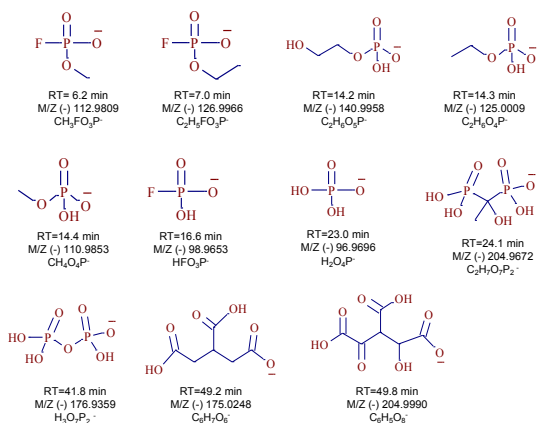
Figure 7. MS Base Peak Chromatogram of Sample 3



**FIGURE 7. Components Identified from Sample 3**

| Components Identified in Cycle Aged Exhibited 20% Loss in Capacity Sample 3 |          |          |             |           |  |
|---|----------|----------|-------------|-----------|--|
| Peak #  | RT (min) | m/z      | Formula (-) | Delta ppm | Name (Based on MS results)*                        |
| 1   | 3.2-3.6  | 125.0009 | C2H6O4P     | 0         | Phosphate Esters                                   |
|   |          | 155.0116 | C3H8O5P     | 0.6       |  |
|   |          | 169.0272 | C4H10O6P    | 0.6       |  |
|   |          | 185.0222 | C4H10O6P    | 0.6       |  |
| 2   | 3.8      | 139.0166 | C3H8O4P     | 0.4       | Phosphoric acid                                    |
| 3   | 4.0      | 89.0244  | C3H5O3      | 0.1       |  |
| 4   | 4.2      | 75.0088  | C2H3O3      | -0.2      | Methyl Carbonate                                   |
| 5   | 5.3      | 139.0071 | C3H7O4S     | 0.4       | Propyl sulfate                                     |
| 6   | 6.2      | 112.9810 | CH3O3FP     | 0.3       | Methyl phosphorofluoridate                         |
| 7   | 7.1      | 126.9966 | C2H5O3FP    | 0.1       | Ethyl phosphorofluoridate                          |
| 8   | 8.2      | 123.0122 | C3H7O3S     | 0.3       | Propyl sulfonate                                   |
| 9   | 9.8      | 140.9864 | C2H5O5S     | 0.3       | 2-hydroxyethyl sulfate                             |
| 10  | 10.0     | 155.0020 | C3H7O5S     | -0.3      |  |
| 11  | 14.2     | 140.9958 | C2H6O5P     | 0.1       | 2-hydroxyethyl hydrogen phosphate                  |
| 12  | 14.3     | 125.0009 | C2H6O4P     | -0.1      | ethyl hydrogen phosphate                           |
| 13  | 14.4     | 110.9853 | CH4O4P      | -0.2      | methyl hydrogenphosphate                           |
| 14  | 15.3     | 131.0350 | C5H7O4      | 0         | 3-carboxy-2-methylpropanoate                       |
| 15  | 15.6     | 117.0193 | C4H5O4      | 0.2       | methyl malonate                                    |
| 16  | 15.7     | 133.0143 | C4H5O5      | 0.2       | 3-carboxy-3-hydroxypropanoate                      |
| 17  | 15.9     | 117.0194 | C4H5O4      | 0.2       | succinate  |
| 18  | 16.2     | 103.0037 | C3H3O4      | 0.3       | 2-carboxyacetate                                   |
| 19  | 16.6     | 98.9653  | HO3FP       | 0         | hydrogen phosphorofluoridate                       |
| 20  | 17.1     | 118.9986 | C3H3O5      |           |  |
| 21  | 17.2     | 96.9601  | HO4S        | -0.2      | hydrogen sulfate                                   |
| 22  | 23.0     | 96.9696  | H2O4P       | -0.5      | dihydrogen phosphate                               |
| 23  | 24.2     | 204.9674 | C2H7O7P2    | 0.7       | hydrogen (1-hydroxy-1-phosphono-ethyl)-phosphonate |
| 24  | 24.4     | 190.9517 | CH5O7P2     | 1.1       | Methyl trihydrogen diphosphate                     |
| 25  | 26.1     | 131.0349 | C5H7O4      | -0.4      |  |
| 26  | 26.7     | 175.0249 | C6H7O6      | 0.3       |  |
| 27  | 26.9     | 147.0299 | C5H7O5      | 0.3       | 4-carboxy-3-hydroxybutanoate                       |
| 28  | 27.8     | 161.0092 | C5H5O6      | 0.1       | Ethandetricarboxylate                              |
| 29  | 27.9     | 103.0037 | C3H3O4      | 0         |  |
| 30  | 41.7     | 176.9360 | H3O7P2      | 0.3       | Trihydrogen diphosphate                            |
| 31  | 49.2     | 175.0249 | C6H7O6      | 0.3       | Tricarballic acid                                  |
| 32  | 49.8     | 204.9312 | C6H5O8      | 0.1       |  |

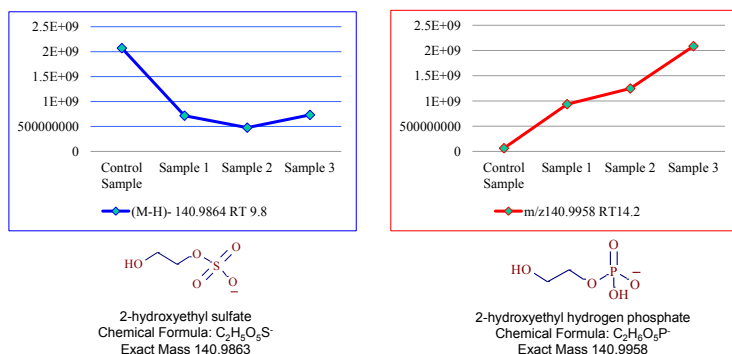
**FIGURE 7. Components Identified from Sample 3**



## Components Profile and Battery Performance

Figure 7 shows the intensity variations in different anode samples of two compounds selected. Because these four anodes were aged under different conditions, the correlation between the compounds and their intensities could be used as an indicator for battery performance.

**FIGURE 7. The Intensity Variations of Two Selected Compounds**



## Conclusion

- Ion chromatography coupled with a Q Exactive mass spectrometer provides a powerful platform for li-ion battery anode impurity and degradant analysis.
- Phosphate degradation products in three aged lithium ion batteries were identified.
- This IC-HRMS platform can be used for QA/QC for lithium-ion battery manufacturers and performance evaluations.
- Further study will be conducted to investigate the relationship of the compounds identified and their intensities with the batteries performance.

## Reference

- [1] The LIB anode samples were provided by a major transportation company.
- [2] See Rosanne Slingsby's AABC2015 poster for in depth IC analysis.

## Acknowledgements

The authors would like to thank the transportation company for providing the lithium ion battery anode samples.

[www.thermofisher.com](http://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  
**Australia** +61 3 9757 4300  
**Austria** +43 810 282 206  
**Belgium** +32 53 73 42 41  
**Brazil** +55 11 3731 5140  
**Canada** +1 800 530 8447  
**China** 800 810 5118 (free call domestic)  
 400 650 5118

**Denmark** +45 70 23 62 60  
**Europe-Other** +43 1 333 50 34 0  
**Finland** +358 10 3292 200  
**France** +33 1 60 92 48 00  
**Germany** +49 6103 408 1014  
**India** +91 22 6742 9494  
**Italy** +39 02 950 591

**Japan** +81 6 6885 1213  
**Korea** +82 2 3420 8600  
**Latin America** +1 561 688 8700  
**Middle East** +43 1 333 50 34 0  
**Netherlands** +31 76 579 55 55  
**New Zealand** +64 9 980 6700  
**Norway** +46 8 556 468 00

**Russia/CIS** +43 1 333 50 34 0  
**Singapore** +65 6289 1190  
**Sweden** +46 8 556 468 00  
**Switzerland** +41 61 716 77 00  
**Taiwan** +886 2 8751 6655  
**UK/Ireland** +44 1442 233555  
**USA** +1 800 532 4752  
 PN64526-EN 08/16S

**Thermo**  
 SCIENTIFIC

A Thermo Fisher Scientific Brand