

Content Discharge



•Detail the Li-ion Battery industry drivers & trends

•Our position in industry and our interest in the application

- Battery research overview
- •How the LiB works and targeted research problems
- Application capabilities
- •Example LiB solutions to tough problems
- •Sample Discussions/LiB Inquiries



Global Drivers

Drivers

Safety

- Consumers Needs
- •Emissions control
- •Globalization in Regulatory
- •Funding, investments, infrastructure
- •Growing performance needs

Result

•Reliable Sustainable energy

- Controllable technology
- •Predictable energy storage and generation

Global Market Objectives

- •Reduce environmental hazard
- •Reduce cost to manufacture
- Increase power & energy of LiB
- Breakdown monopoly
- •Enhance warranties and consumer confidence

Market Influence

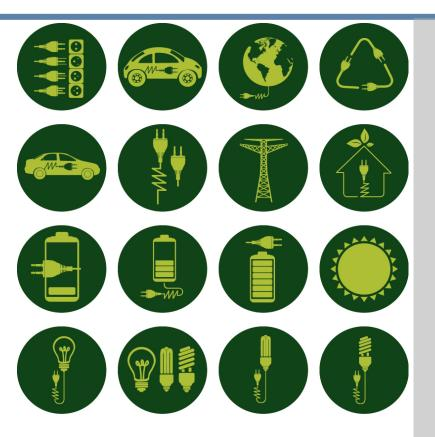
Pervasive technologies- Auto
Plateau of available LiB power
Limit to fuels and alternative energy resources
Start-ups, JVs, regional shifts

Forecast

End market CAGR 23.3% 2013-2020
By 2013 >100 competitive players in LiB
EV market to be 2X consumer LiB by 2023
Strong funding of academic, contract labs, and government funded test-houses



Our Responsibility Behind Understanding LiB



We have the responsibility to add our experience in the areas of research that make our world healthier, cleaner, and safer.

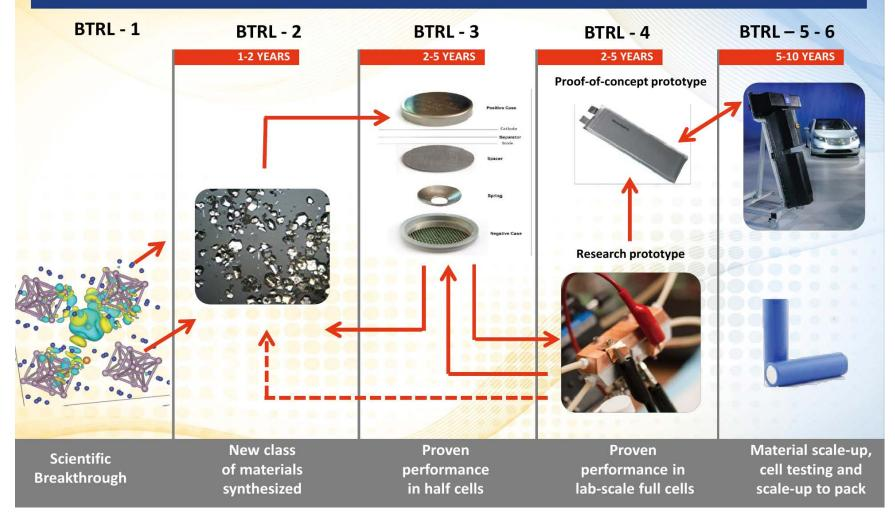
Build new Solutions that take a closer look at the materials, changes, and future of battery analysis.

Help our global partners achieve their long term research goals!



The Ten Year Cycle of R&D to Market

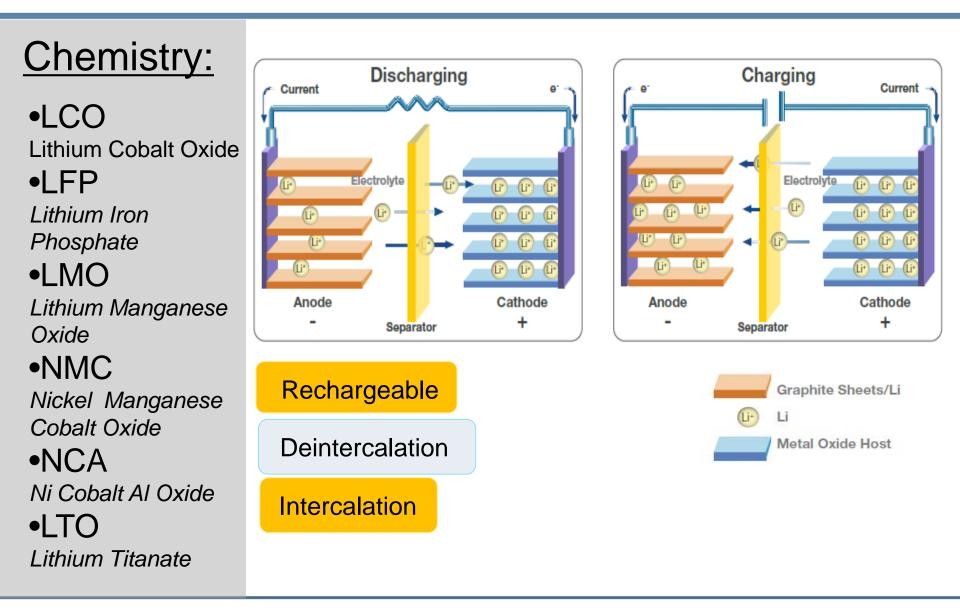
Battery Technology Readiness Level (BTRL)





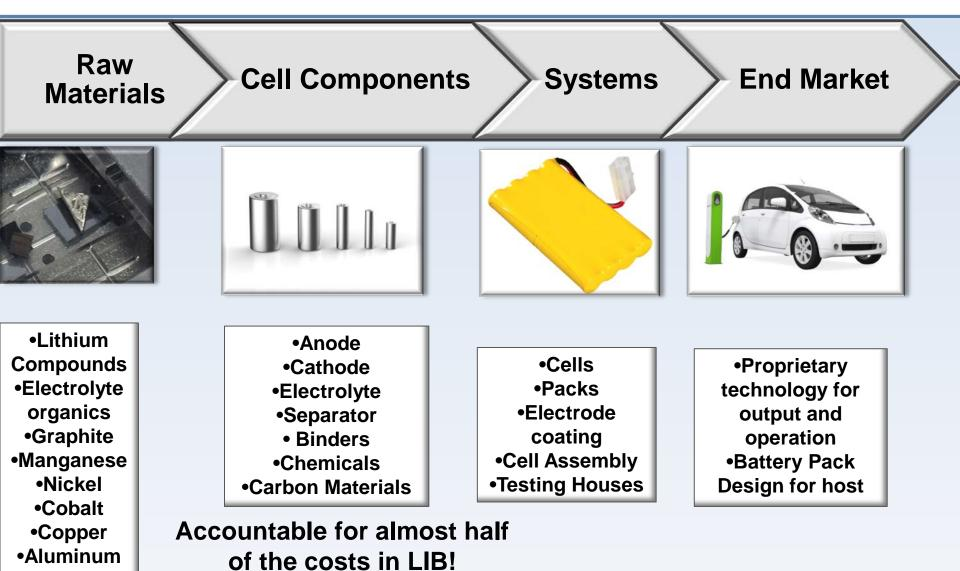
Property of Argonne National Laboratory: https://anl.app.box.com/batterytechnologyreadiness

How the Lithium-ion Battery Works





LiB Value Chain

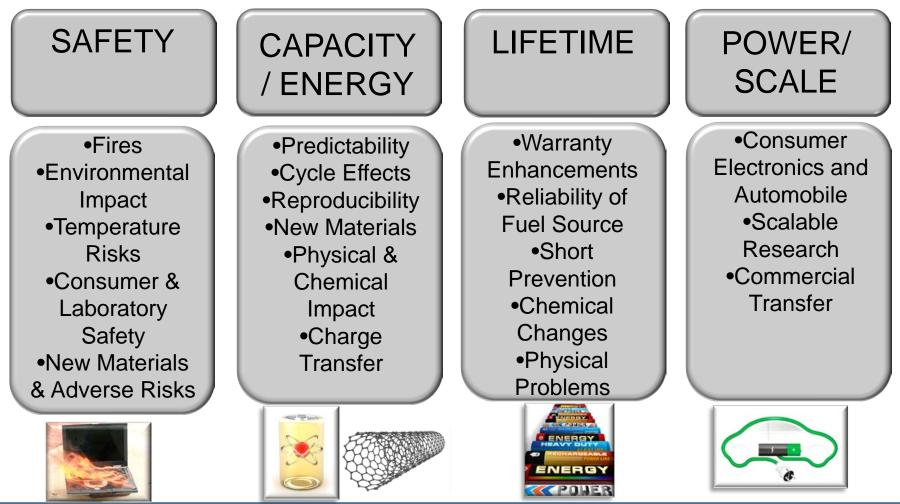


Thermo Fisher

SCIENTIFIC

Research Tradeoffs

Balance between tradeoffs is critical to successful research! How do we maximize the balance?



Thermo Fisher

Key Problem Areas- The Battery Breakdown

Can Steel Plate Wave Spring Cathode (+) Anode (-) Separator & Electrolyte		Discharge e - Discharge Li+ Li+ Anode (-) Cathode (+) Separator & Electrolyte				
Cell	Electrolyte	Separator	Binder	Anode	Cathode	
 In situ activity Optimum cell chemistry Prevention of leaks Internal Impedance Stability of varying components and cell 	 Additives Ion dispersion Gas generation Flammability Low flashpoint Breakdown products 	 Porosity effects Copolymer characterization Resistance Mechanical strength Impurities Thickness Temperature Limits 	 Homogeneity Surface area control Composition Heat resistance Material variance Impurities Viscosity Adhesion 	 Lithium deposition Dissolution Expansion SEI Layer Silicon Behavior Ion Dispersion Particle Morphology 	•Oxide formations •Volume changes •Film growth •Functional group ID •Dendrites •Impurities •Capacity effects	

Define Battery Question and Problems

ex situ



- Battery disassembly to analyze Individual Components Upon
- Inert Sample change/failure. Transfer
- Reactive Analysis
- Destructive analysis (Battery De-assembled)

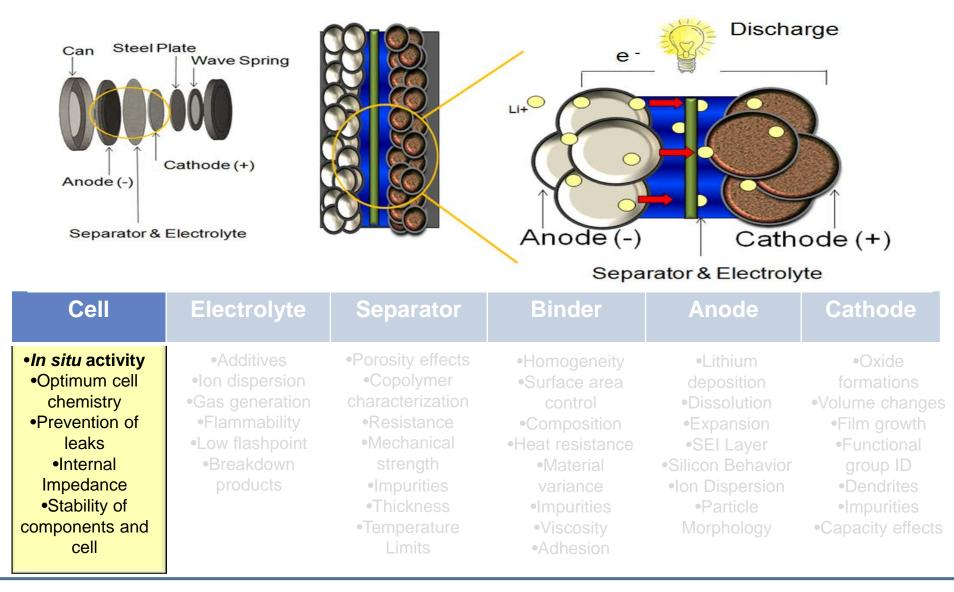




Analysis under load, in a working environment to observe real-time cell activity.

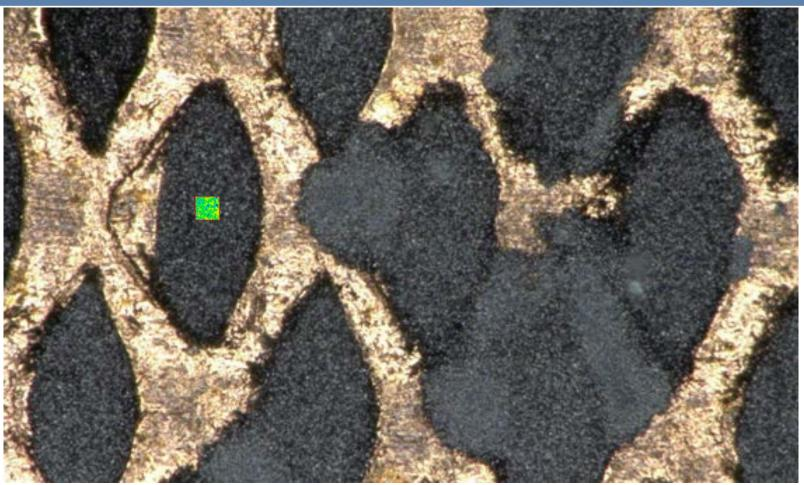
- Can view during operation/charge cycles
- Put into a 'real' situation
- Assemble a battery
- Proactive Analysis

Key Problem Areas- in situ Cell Investigations





in situ Raman: Lithiation of Graphite



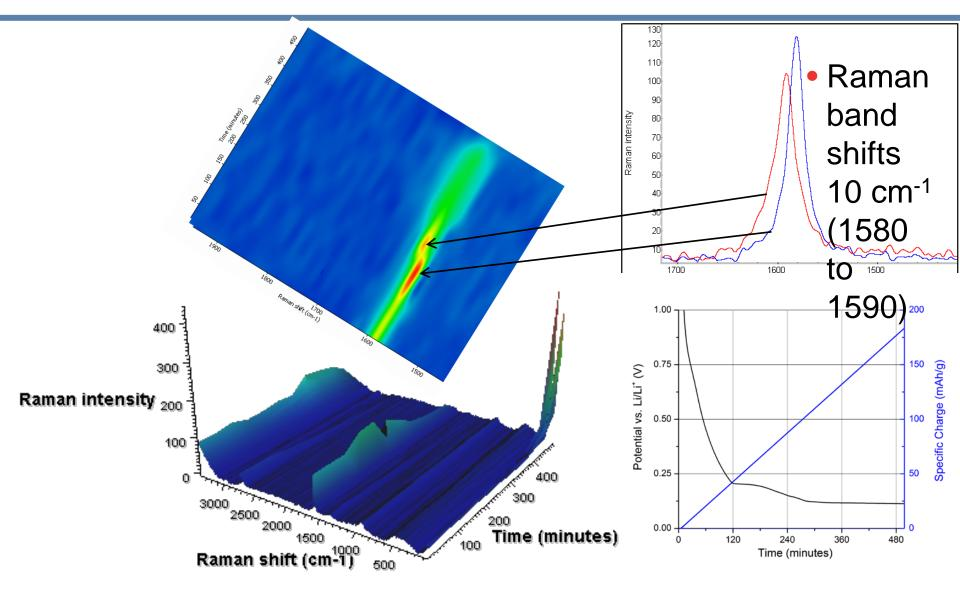
- Graphite coated on wire mesh current collector
- Representative area examined by Raman

Image and electrochemical data provided by EL-CELL, use of ECC-Opto-Std optical electrochemical cell, 2015



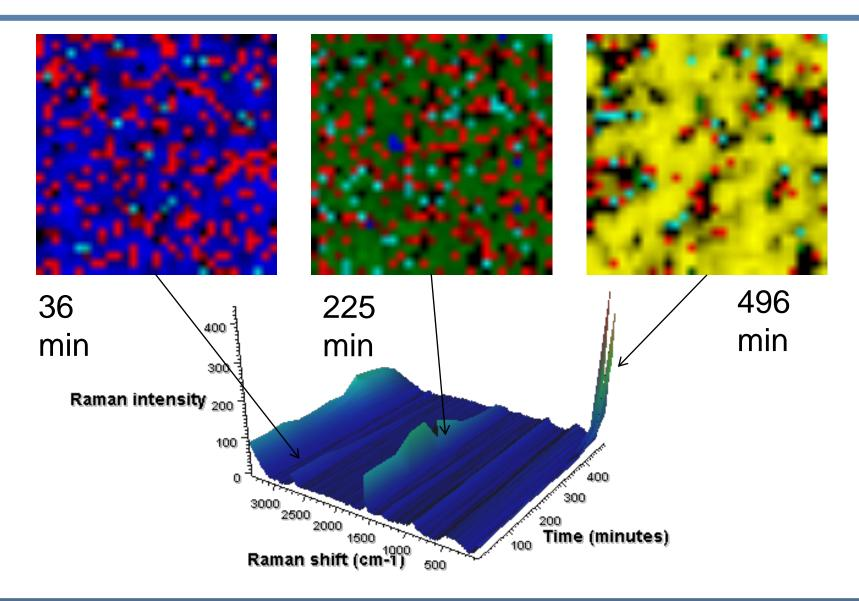


in situ Raman: Change in Spectrum Over Time





in situ Raman: Change in Raman Image Over Time



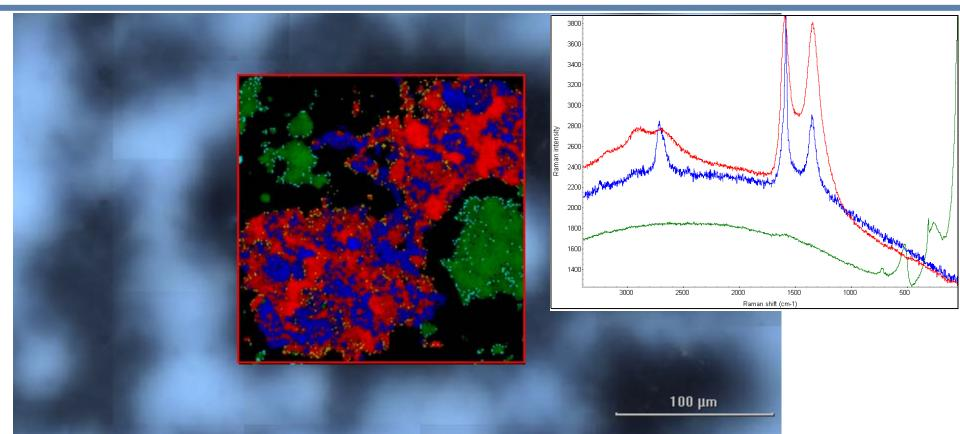


Key Problem Areas- Ex Situ Electrode Investigations

Can Steel Plate Wave Spring () () () () () () () () () () () () ()		e - Discharge Li+ Cathode (+) Anode (-) Cathode (+) Separator & Electrolyte				
Cell	Electrolyte	Separator	Binder	Anode	Cathode	
 In situ activity Optimum cell chemistry Prevention of leaks Internal Impedance Stability of varying components and cell 	 Additives Ion dispersion Gas generation Flammability Low flashpoint Breakdown products 	 Porosity effects Copolymer characterization Resistance Mechanical strength Impurities Thickness Temperature Limits 	 Homogeneity Surface area control Composition Heat resistance Material variance Impurities Viscosity Adhesion 	 Lithium deposition Dissolution Expansion SEI Layer Silicon Behavior Ion Dispersion Particle Morphology 	•Oxide formations •Volume changes •Film growth •Functional group ID •Dendrites •Impurities •Capacity effects	

Thermo Fisher

Post Diagnostic Li-ion Battery Anode, 2yr Cycle

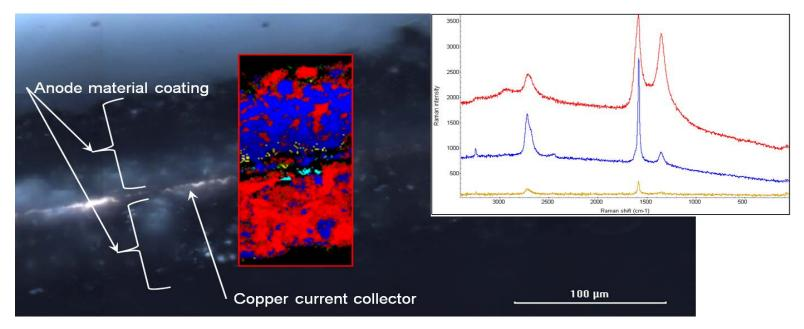


 Electrode material (dark areas) lies between overlayer of separator particles (light areas)

- Red (29% area) & blue (20% area) are variations in SEI layer
- Green is separator particle

Ex situ Analysis of a Cross-Sectioned Anode Material

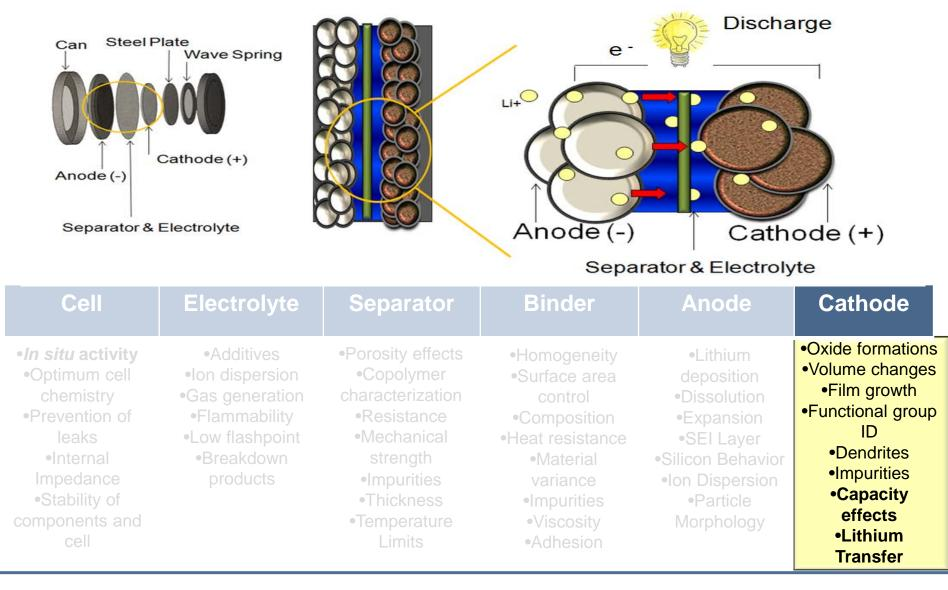
Cross-sectioned anode material in the ex situ transfer cell



The red color indicates the presence of carbon black while the blue color represents graphite. The distribution of these materials on the two sides of the electrode is significantly different. The copper current collector is in the center.

50X long working distance objective, 532 nm laser (2.0 mW), area imaged 76 μ m x 160 μ m, image pixel size 1 μ m, 0.2 s exposure time, 4 scans

Key Problem Areas- Surface analysis of Cathodes



Thermo Fisher

K-Alpha⁺ Sample Transfer Capability and weak signal detection

Analysis Examples:

<u>Electrodes</u>

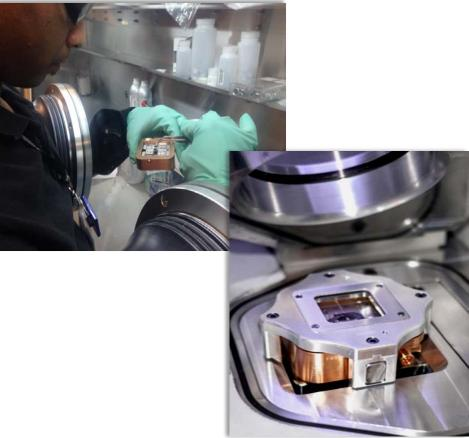
- Surface characterization of pristine material
 - Confirm oxidation state, composition, Li gradients
- Ex-situ characterisation after cycling
 - Composition & variation with depth of SEI
 - Variation in surface composition of electrode material

Separators

- Surface characterization of pristine material
 - Confirm surface chemistry
- Ex-situ characterisation after cycling
 - Look for polymer degradation
 - Deposition of material from electrodes & electrolyte

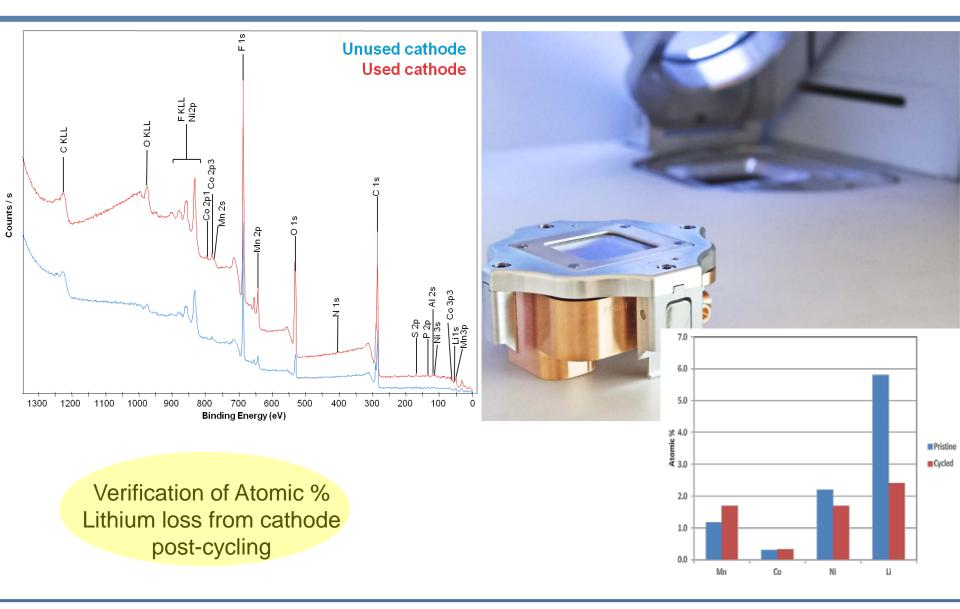
Inert atmosphere transfer

- Load samples in glove box
- Transfer under vacuum to K-Alpha⁺



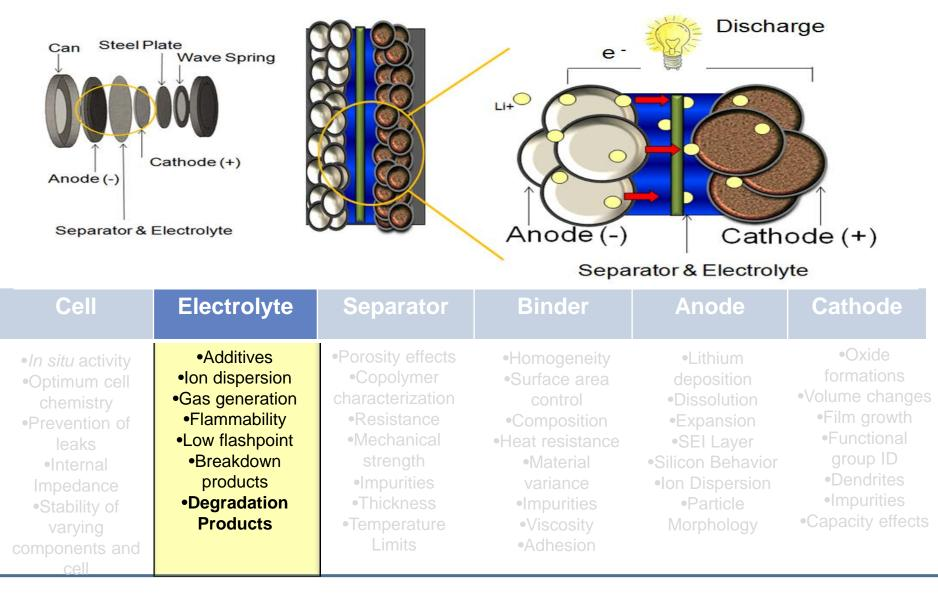


XPS Comparison of Pristine and Cycled Cathode



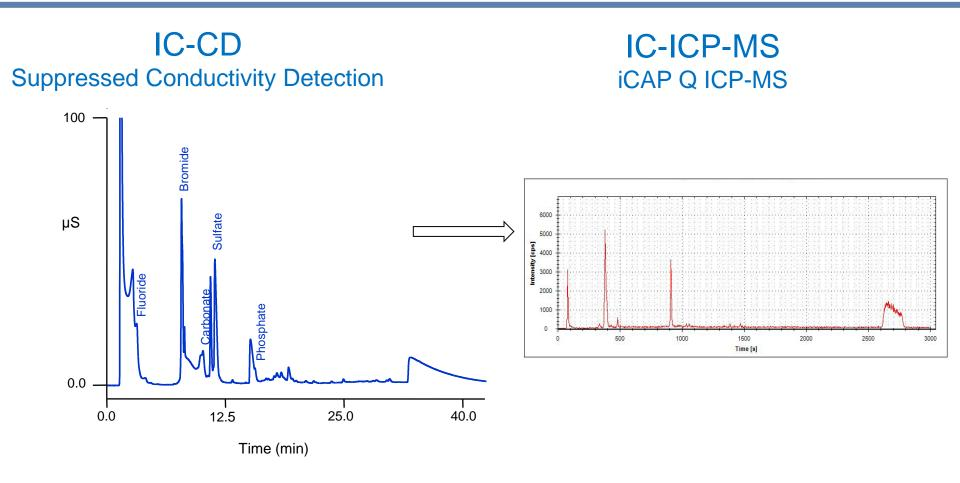


Key Problem Areas- Electrolytes



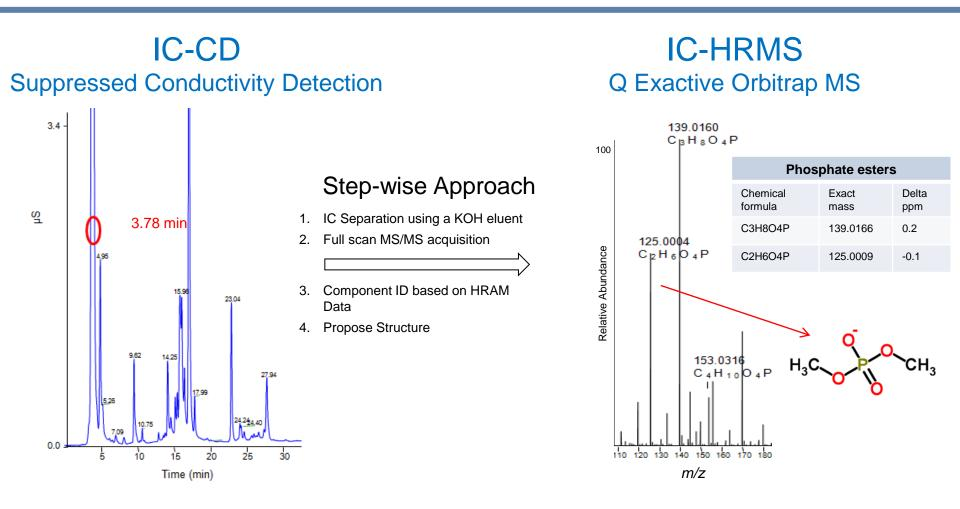
Thermo Fisher SCIENTIFIC

Example 1: Li-Ion Battery Analysis: IC-ICP-MS



Analyze ³¹P Containing Products in the Presence of Other Elements

Example 2: Li-ion Battery Analysis: IC-HRMS

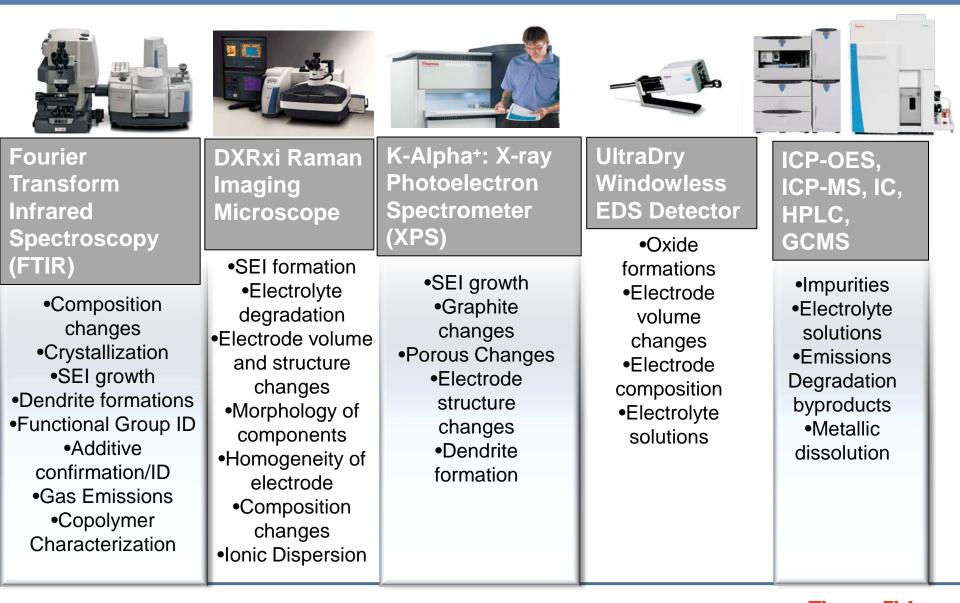


Component Identification in Untargeted and Unknown Workflows

Source for Dimethyl phosphate image: CSID:2982799, http://www.chemspider.com/Chemical-Structure.2982799.html (accessed 00:59, Feb 5, 2015)



Thermo Scientific Building Block of Application Solutions

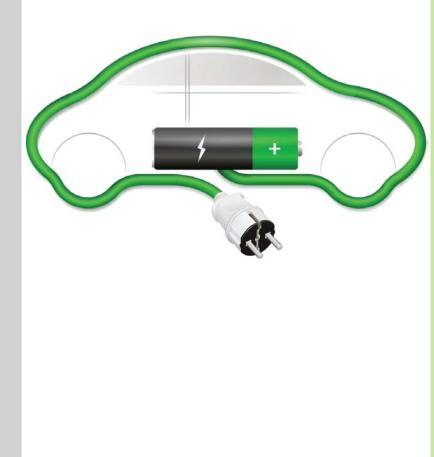


Thermo Fisher

The Automotive and Advanced Materials Industry

Targeted Pieces

- •Carbon Materials
- •Glass technology
- •Display Materials
- •Lighting Advancements
- •PowerTrain
- •New Energy
- •Renewable Energy
- •Secondary Batteries (Li-ion)
- •Carbon Filled Materials
- •Brake Advancements
- Lubricants
- Rubbers
- Adhesives
- Recyclables
- Plastics
- •Paints & Coatings
- •Laminates



Goals in Research

- •Materials: Weight
- Reduction
- •Materials: Strength
- Enhancement
- Catalysis
- •Scratch Resistance
- •Safety
- •Failure Analysis
- •Fire Prevention (Flame Retardants)
- Emissions Control
- •Lifetime of Product
- •Corrosion
- Resistance
- Corrosion Resistance
- •Reduce, Reuse,
- Recycle
- Heat Resistance
- Color Retention

