

High-quality *in operando* X-ray diffraction analysis of pouch bag lithium-ion batteries



Introduction

Lithium-ion batteries have become increasingly important in key technologies of our day-to-day lives. Especially the long-term performance of batteries is crucial for their applications in new markets such as energy storage in electric vehicles or even large-scale electric storage devices combined with photovoltaic energy sources.

For these large-scale applications a key requirement is very good long-term and cycling performance. Improvement of this performance requires studying ongoing aging processes inside batteries continuously and without the need to destroy and disassemble them prior to the analysis (*in operando* analysis). This is especially important as many cell components deteriorate due to the disassembly process or when exposed to air or moisture.

Often X-ray synchrotron or neutron radiation based methods are used in this context [1-4]. The strengths of these methods are their high penetration capabilities and their ability to use specially encapsulated samples or even industrially manufactured samples directly. Nevertheless, a drawback of these methods is the restricted access as they are only available at large-scale research facilities and obtaining the required beam time is typically a longer process.

While neutron scattering is very sensitive to light elements such as Li, X-ray powder diffraction provides a much higher resolution of the cell parameters and therefore both are complementary techniques. A large variety of experimental cells for XRD studies has been developed such as specially adapted coin cell type batteries for reflection geometry. Typically XRD setups of lab diffractometers are based on using copper radiation and rely on using so-called 'half-cells', containing only one active electrode while the other electrode is then only a thin foil of lithium metal.

Although this is adequate to study the principal transformation mechanism and the corresponding structural changes of the electrode active material via diffraction methods, it is insufficient for studying complex aging processes. These depend heavily on even the smallest changes in cell chemistry and material composition and require monitoring the structural evolutions of both electrodes at the same time. For such aging studies one cannot rely on half-cell studies but has to observe the cells under real operating conditions in their native environment.

The pouch bag cell design offers this possibility. It is easy to manufacture reproducibly. It is versatile enough for experimental changes and offers the possibility to use all standard battery grade materials and components. An industrial prototype can be assembled

Summary

Pouch cell batteries can be used in combination with hard X-rays in transmission geometry to perform *in situ* and *in operando* studies during charge/discharge cycles and for aging studies. It is then possible to correlate variations in the crystallographic structure of the elements in the cell directly with the amount of Li incorporated in the electrodes. Moreover, the efficiency of the GaliPIX^{3D} detector can considerably improve data quality compared to standard Si-based detectors.

for the continuous monitoring of cells subjected to different aging procedures. In synchrotron beam lines such pouch cells have already been used successfully [5].

The battery research group headed by Dr. habil. Ralph Gilles (TU München) at the Heinz Maier-Leibnitz Zentrum (MLZ) in Garching (Germany) has adapted and further developed this method in order to use standardized pouch cells for diffraction experiments on a lab XRD instrument like the Emyrean using hard X-ray radiation. This allows investigation of the structural dynamics of fully closed pouch bag type batteries as used as industrial prototype samples during typical aging procedures directly and without any prior dismantling.

The data and method described in this application note resulted from a common work of Dr. Stefan Seidlmayer (TU München) and Armin Kriele (Helmholtz Zentrum Geesthacht) from the Heinz Maier-Leibnitz Zentrum, MLZ in Garching (Germany).

The MLZ is a neutron research facility open for user experiments in science and industrial application

Measurement setup

The first set of measurements was performed at the Heinz Maier-Leibnitz Zentrum (MLZ) in Garching (Germany), where a customized transmission setup was used for the *in situ* measurements, using a PANalytical Empyrean diffractometer with an X'Celerator detector.

The second set of measurements was performed at PANalytical headquarters in Almelo (the Netherlands), and it includes a comparison of the results obtained with an X'Celerator and the GaliPIX^{3D} detector.

Mo radiation was used for both measurements, in combination with a focusing beam X-ray mirror. Photographs of the setups used and the pouch cell are presented in Figure 1.

The cell was cycled twice between discharged and charged states following a sequential CCCV (constant current – constant voltage) procedure. First the cell is charged incrementally in ten steps with a C/10 rate and 2 h stops in the current flow. Then the cell is charged further until it reaches a certain cut-off voltage potential and the current flow drops below the cut-off threshold. The cell potential and current flow during the procedure are shown in Figure 2. The *in situ* XRD measurements were performed continuously over several days, and consisted of two complete charge/discharge cycles of the battery cell (see Figure 3). Powder diffraction scans were collected continuously; each scan in the range 5° to 55° 2theta with a measurement time for each scan of 16 minutes 44 seconds. Subsequently, scans measured during the ramping of the current were discarded, and the scans performed at constant zero current value were analyzed. A script was written in HighScore Plus (version 4.1 [6]) to automatically merge all the scans measured at the same conditions.

Instrument setup

Samples were measured with the following parameters:

Instrument	Diffractometer
Radiation	Mo K α
Incident beam optics	Focusing mirror with ¼ Ta divergence slit, ¼ Ta anti-scatter slit, 0.02 rad Soller slit, 10 mm mask
Sample stage	3-axes cradle in combination with customer-designed sample holder
Diffracted beam optics	Anti-scatter slit, 0.02 rad Soller slits
Detector	X'Celerator and GaliPIX ^{3D}
Scan parameters	5° - 55° 2 θ scans, 16 min scan time

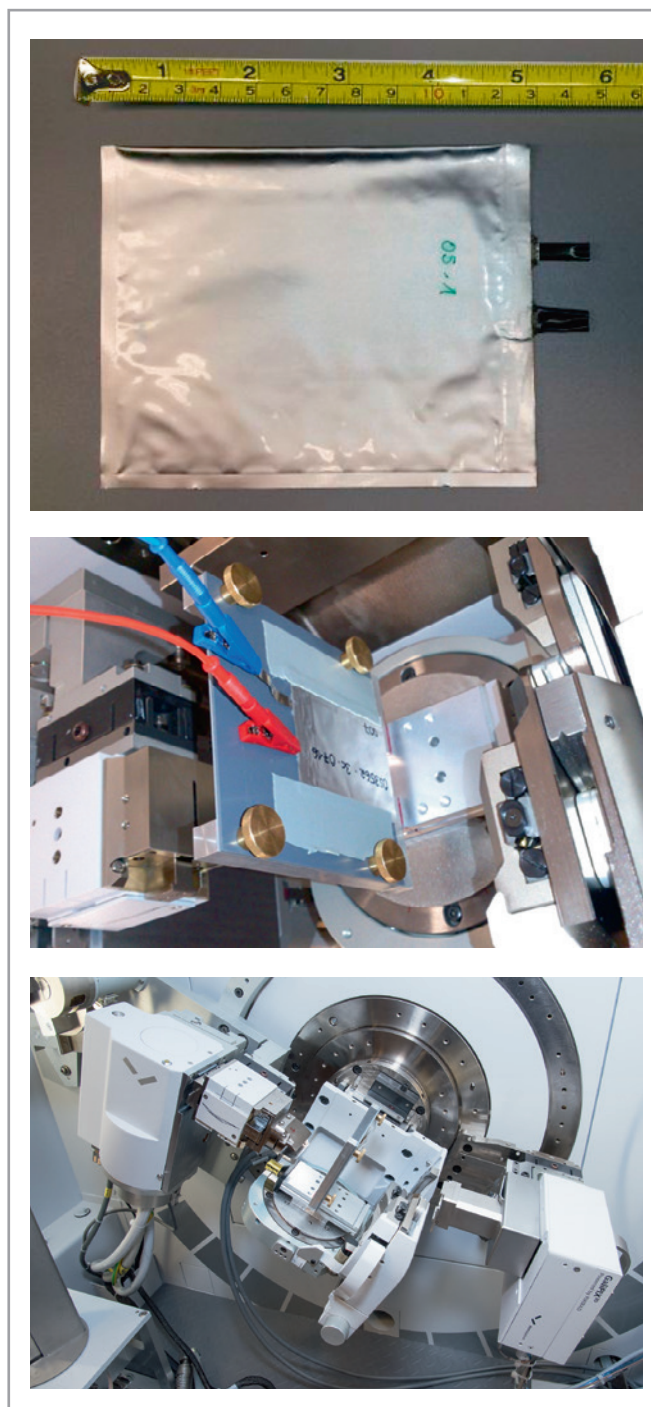


Figure 1. Top to bottom: typical pouch bag battery after assembly; setup used at the Heinz Maier-Leibnitz Zentrum (MLZ) in Garching for *in situ* measurements; measurement setup used for testing GaliPIX^{3D} performance

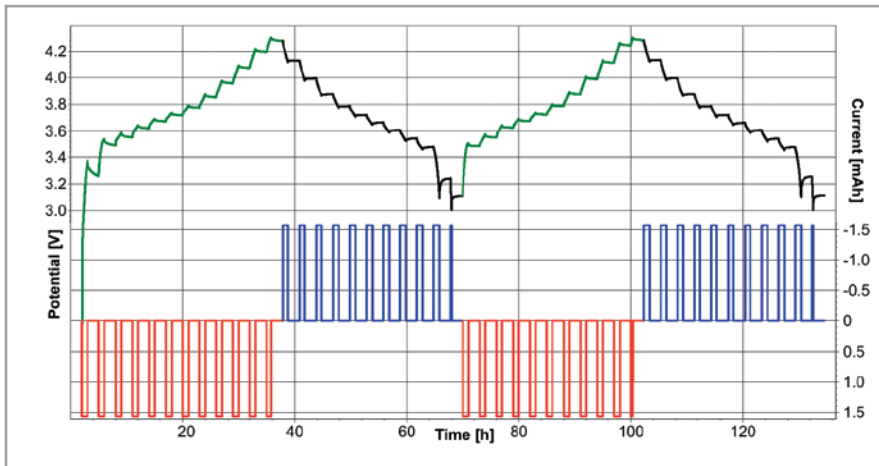


Figure 2. Cell potential and current during XRD measurements. In green/black the potential during charge and discharge respectively (left axis). In red/blue the current (right axis)

Basic principle of a Li-ion battery and pouch bag battery cells

The energy storage in a Li-ion battery is achieved by forcing Li ions to move from the cathode active material to the anode active material by applying an electrochemical force (Figure 3).

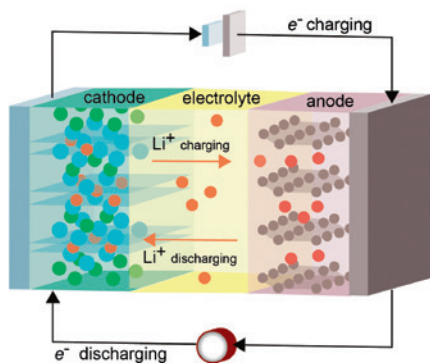


Figure 3. Schematic view of a lithium-ion battery

During charging an external electrical power source applies an over-voltage (a higher voltage than the battery produces, of the same polarity), forcing the current to pass in the reverse direction. The lithium ions then migrate from the cathode to the anode, where they become embedded in the layered electrode material in a process known as intercalation.

Lithium batteries used in the measurements were assembled in the so-called pouch bag format. In this assembly type, the electrodes of the battery are protected inside a thin polymer composite hull (pouch).

This pouch is a complex polymer/metal assembly of an aluminium foil sandwiched between polymer layers. The overall thickness of such a pouch bag is ~100 μm . Figure 4 shows a schematic view of such a cell.

The electrodes were prepared by coating thin metal current collector foils with dispersions of active materials in NMP (N-methyl-2-pyrrolidone). The dry thickness of the coatings was ~60 μm .

For anodes, a 10 μm Cu foil was coated with artificial graphite and PVDF-binder.

For cathodes, a 15 μm Al foil was coated with $\text{LiNi}_{0.33}\text{Mn}_{0.33}\text{Co}_{0.33}\text{O}_2$ (LiNMC) active material mixed with PVDF-binder and carbon black as a conductive additive. The electrodes are separated by a 25 μm thin layer of electrically insulating polypropylene material.

After stacking the electrode layers in the pouch bag a standard electrolyte solution of LiPF_6 dissolved in a mixture of EC:EMC (ethylene carbonate: ethyl methyl carbonate) was added prior to heat-sealing the pouch bag cell.

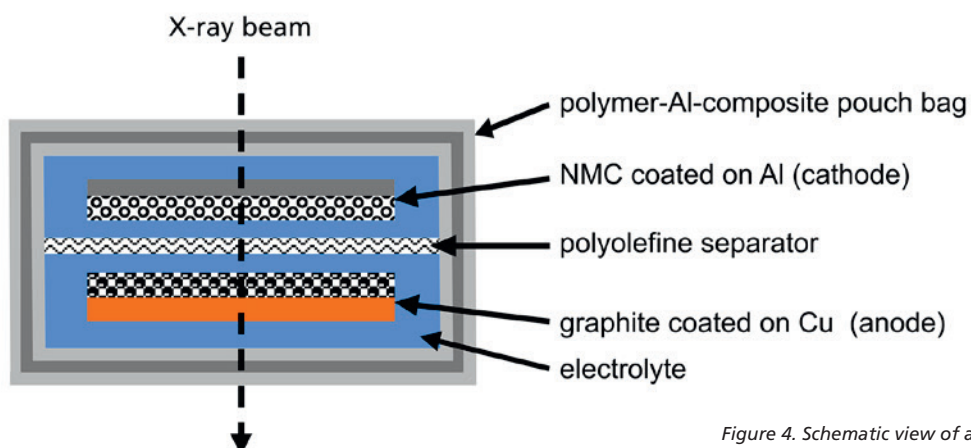


Figure 4. Schematic view of a pouch bag cell

Results and discussion

Automatic Rietveld and Pawley mixed fits were performed with HighScore Plus on every scan, to obtain the cell parameters for the LiNMC and the graphite phases. An example of the fit is shown in Figure 6.

The cell parameters obtained during the fit were automatically exported to a spreadsheet file, and the results can be found in Figure 5.

From the automated fits changes in the lattice parameters can be observed, which relate directly to the structural evolution of the electrode materials during the intercalation/deintercalation process. Most striking feature is the different behavior of the a - and c -axis of the LiNMC cathode. During the charge process, which corresponds to removal of lithium ions from the structure of the cathode, the a cell parameter first shows a decrease followed by a shorter region where it is nearly constant. On the other hand the c -axis shows first a strong increase and then switches to a decrease just before the cell is fully charged. Both these trends are reversible upon discharge of the battery.

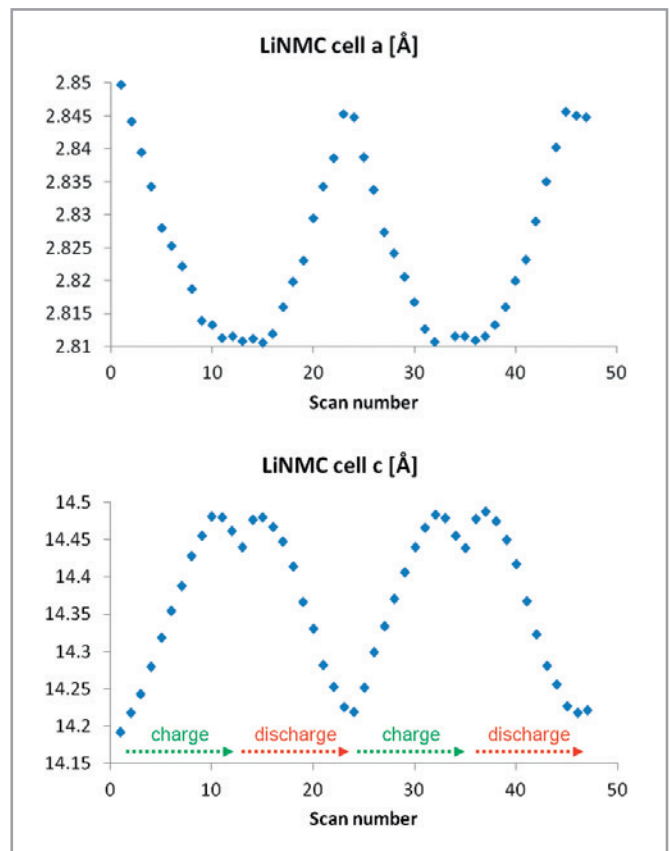


Figure 5. Fitted LiNMC lattice parameters

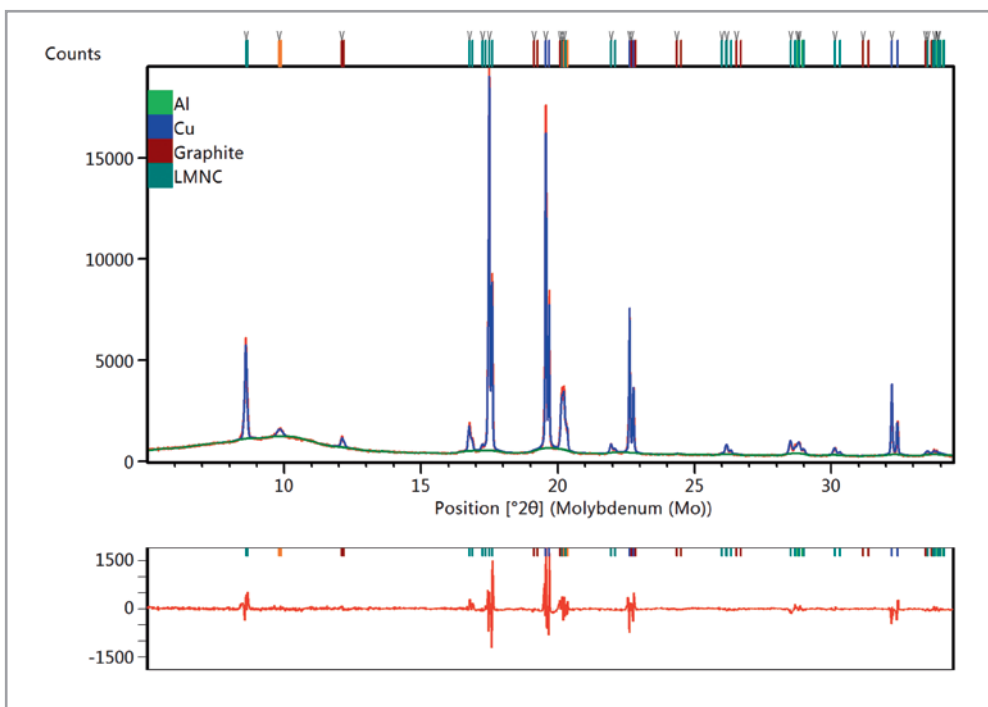


Figure 6. Example of the data including the different phases, the automatic fit and the difference plot in the lower panel

Results and discussion (continued)

The structure of the layered transition metal oxide LiNMC is shown in Figure 7.

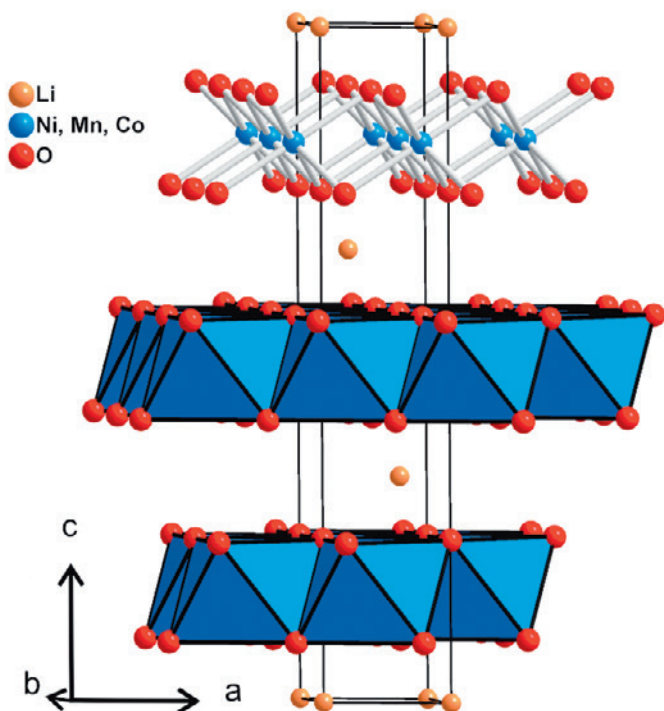


Figure 7. Projection of the crystal structure of the LiNMC cathode along b

The transition metal atoms are arranged in layers of edge-sharing MO_6 octahedra. These layers are stacked along the c -axis with an ABC-packing sequence. The lithium atoms are inserted in the interlayer space between the MO_6 -layers and are octahedrally coordinated by oxygen atoms. These observations can be rationalized considering the structural features. During charging the lithium atoms leave the structure and an increase in the oxygen-oxygen electrostatic repulsion expands the lattice along the c -axis. At the same time the average oxidation state of the transition metal atoms is successively increasing while Li^+ ions are leaving the structure to maintain the overall charge balance. This leads to a stronger ionic interaction between metal and oxygen atoms leading to a shortening of the M-O bonds which are lying within the a - b -plane and thus resulting in a decrease of the a lattice parameter.

When the sample mass and thus the absolute number of moles of active LiNMC material in the cell is known, one can also calculate the amount of $x(Li)$ in the formula as every transferred electron equals one lithium atom. This means that for the charge/discharge process one can also write the formula of LiNMC as $Li_{1-x}Ni_{0.33}Mn_{0.33}Co_{0.33}O_2$. It is interesting to plot either the lattice parameters or the parameter ratio c/a versus the calculated amount of $x(Li)$ in $Li_{1-x}NMC$ as shown in Figure 8.

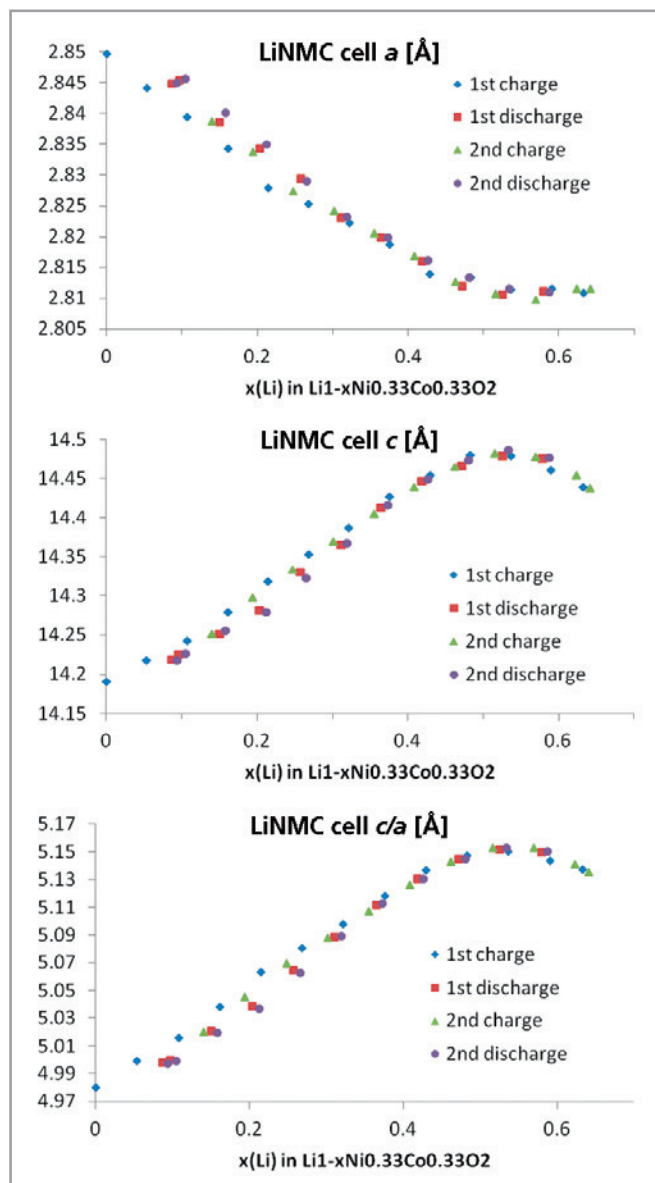


Figure 8. Cell parameters obtained for the LiNMC and graphite phases, as a function of the Li content

Results and discussion (continued)

This correlation shows that there is some hysteresis between the first and successive charge/discharge cycles in regard to the lattice parameters. This effect is called irreversible capacity loss (ICL). It is a result of the very slow uptake of lithium ions upon re-insertion into the host $\text{Li}_{1-x}\text{NMC}$ structure when reaching x -values close to 1. By measuring the linear correlation between the lattice parameter ratio c/a and the lithium content $x(\text{Li})$ during the 1st and 2nd cycle of the cell with *in situ* XRD it is also possible to obtain a value for the

amount of active lithium in the cathode. It is important to highlight that this value can then be compared to the electrically observable capacity. Deviations of these two values from each other give valuable insights for the understanding of aging processes in lithium-ion batteries [7].

By using the isolines plot available in HighScore, it is also possible to observe the variation of the cell parameters directly by following the shifting of the Bragg peaks for LiNMC as well as for graphite, see Figure 9.

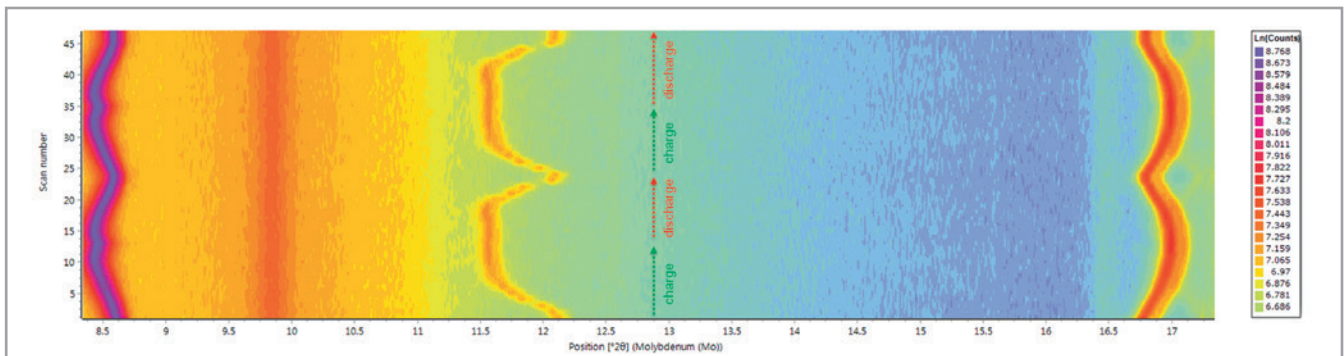


Figure 9. Variation in the XRPD data during two full charge and discharge cycles. The peaks at about 8.5° and 17° 2theta belong to the LiNMC phase, (003) and (101) reflections respectively. The peak at about 12° 2theta is the (002) reflection of graphite.

With the transmission geometry setup and the pouch bag cells in the *in situ* XRD setup, it is not only possible to follow the structural changes due to the deintercalation of lithium from LiNMC on the cathode. Simultaneously the changes due to the lithium intercalation into the graphite host lattice on the anode side of the cells can be followed (see Figure 10).

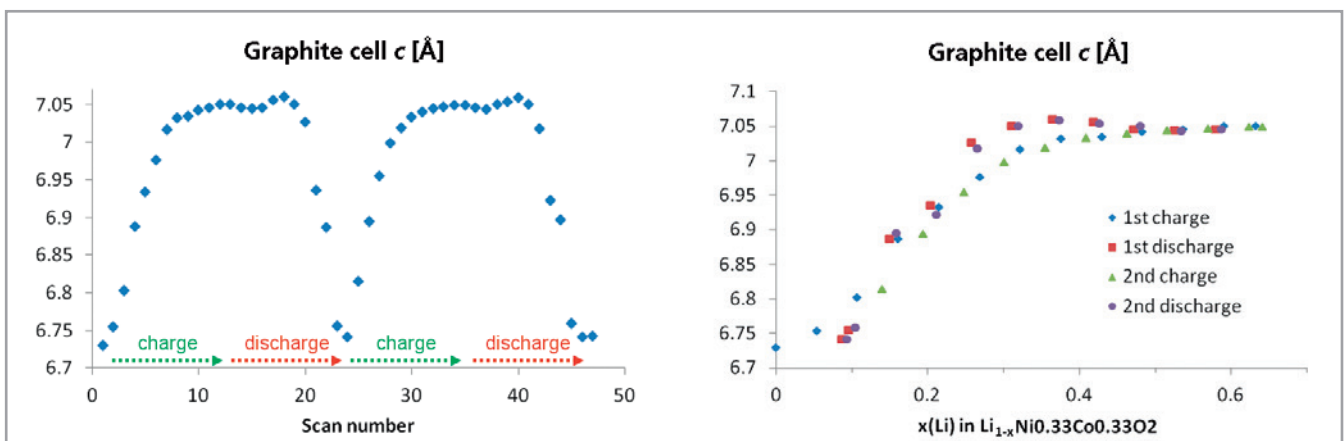


Figure 10. Graphite cell parameters as a function of scan number (left) and lithium content (right)

The detailed structural dynamics involved in the transformation of the structure of pure graphite into the structure of the fully intercalated graphite, LiC_6 , can also be observed. The graphite structure consists of graphene layers packed along the c -axis.

Following the lattice parameters (Figure 10) starting with a graphite basic structure, first a lattice expansion along the c -axis can be observed. This is explained with the insertion of the lithium atoms into the interlayer space of the graphite host structure resulting in an expansion along the c -axis to accommodate the additional lithium ions in the structure.

Improving the data further

The GaliPIX^{3D}, a 1D/2D detector with a high-quality CdTe sensor, is the ideal detector for this application. Over the X'Celerator used in this study it offers a ~five times higher intensity which can be used to collect data with higher resolution (by narrowing slits) or in the shortest possible time frame. The GaliPIX^{3D} can also be used for the analysis of amorphous or highly disordered materials using the pair distribution function (PDF) method and/or transmission X-ray imaging and computed tomography (CT).

Figure 11 shows the performance difference between the two detectors on the same pouch cell. It clearly demonstrates the benefit of using GaliPIX^{3D} for *in operando* measurements on complete electrochemical cells in the transmission geometry.

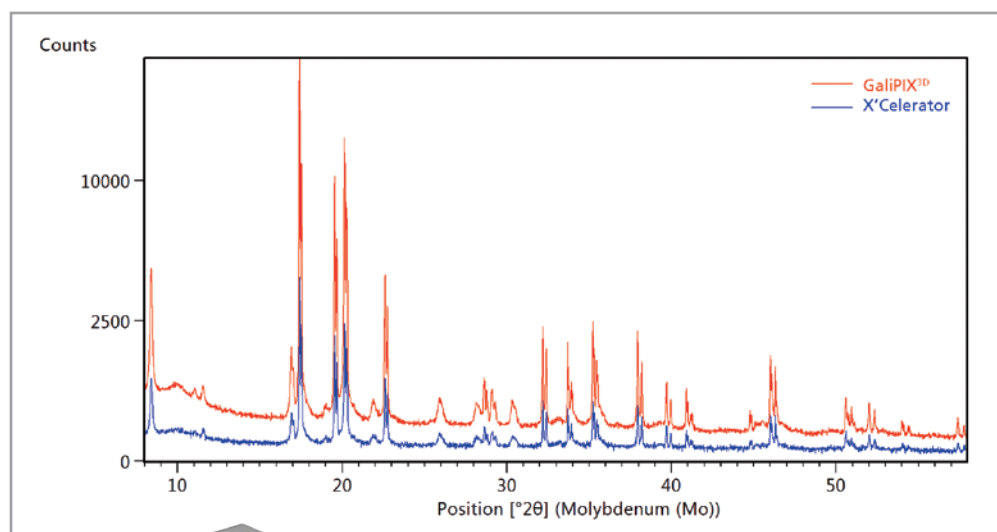


Figure 11. Comparison of data measured using the same setup and total scan time with GaliPIX^{3D} data (red) and X'Celerator data (blue)



- [1] V. Zinth, S. Seidlmayer, N. Zanon, G. Crugnola, M. Schulz, R. Gilles, and M. Hofmann, *Journal of The Electrochemical Society*, **162** (3), A384-A391 (2015).
- [2] M. Morcrette, Y. Chabre, G. Vaughan, G. Amatucci, J. B. Leriche, S. Patoux, C. Masquelier, and J. M. Tarascon, *Electrochimica Acta*, **47** (19), 3137-3149 (2002).
- [3] S. Seidlmayer, J. Hattendorff, I. Buchberger, L. Karge, H. A. Gasteiger, and R. Gilles, *Journal of The Electrochemical Society*, **162** (2), A3116-A3125 (2015).
- [4] C. Kortenbruck, R. Gilles, A. Senyshyn, V. Zinth, M. Hofmann, C. von Lüders, O. Dolotko, M. Mühlbauer, S. Seidlmayer, and N. Paul, *Neutron News*, **26** (1), 29-32 (2015).
- [5] K. W. Nam, W. S. Yoon, H. Shin, K. Y. Chung, S. Choi, and X. Q. Yang, *Journal of Power Sources*, **192** (2), 652-659 (2009).
- [6] T. Degen, M. Sadki, E. Bron, U. König and G. Nénert. *Powder Diffraction*, **29**, S13-S18, (2014)
- [7] I. Buchberger, S. Seidlmayer, A. Pokharel, M. Piana, J. Hattendorff, P. Kudejova, R. Gilles, and H. A. Gasteiger, *Journal of The Electrochemical Society*, **162** (14), A2737-A2746 (2015).

Conclusions

Pouch cells batteries can be measured *in situ* and *in operando* with hard X-rays in transmission geometry as a stand-alone method or as a complementary technique to neutron diffraction. Information about the cell parameters of anode and cathode have been obtained using the HighScore Plus software. GaliPIX^{3D} is the ideal detector for this kind of study, thanks to its extremely high efficiency in detecting hard X-rays.

About PANalytical

PANalytical's mission is to enable people to get valuable insight into their materials and processes. Our customers can be found in virtually every industry segment, from building materials to pharmaceuticals and from metals and mining to nanomaterials. The combination of our software and instrumentation, based on X-ray diffraction (XRD), X-ray fluorescence (XRF) and near-infrared (NIR) spectroscopy as well as pulsed fast thermal neutron activation (PFTNA), provides our customers with highly reliable and robust elemental and structural information on their materials and is applied in scientific research and industrial process and quality control.

PANalytical employs over 1,000 people worldwide. The company's headquarters are in Almelo, the Netherlands. Fully equipped application laboratories are established in Japan, China, the US, Brazil, and the Netherlands. PANalytical's research activities are based in Almelo (NL) and on the campus of the University of Sussex in Brighton (UK). Supply and competence centers are located on two sites in the Netherlands: Almelo (X-ray instruments) and Eindhoven (X-ray tubes), in Nottingham, UK (XRF applications and standards), in Quebec, Canada (fusion sample preparation) and in Boulder CO, US (near-infrared instruments).

PANalytical is active in all but a few countries of the world. This worldwide sales and service network ensures unrivalled levels of customer support.

The company is certified in accordance with ISO 9001 and ISO 14001.

Visit www.panalytical.com for more information about our activities.

PANalytical is part of Spectris plc, the productivity-enhancing instrumentation and controls company.

Access to expertise

With the largest service network we are able to offer the most comprehensive support package possible.

Expertise:

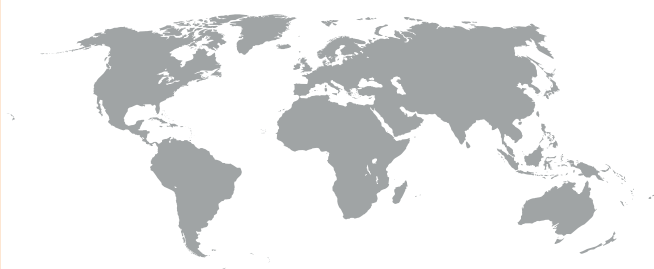
- On-site training available
- XRF training courses
- Performance optimization
- Customizable expertise programs
- Assistance with multi-laboratory standardization

Care Agreements

Our customer support solutions have been developed with your business in mind. They are formulated as a family of four Care Agreements which can be tailored to your specific needs and provide fast, secure and reliable support.

- **ECONOMY:** indispensable coverage for self-sufficient operations
- **ADVANCED:** cost-effective support for routine usage
- **PREMIUM:** flexible package for high equipment usage
- **ELITE:** most comprehensive package for demanding environments

Global and near



PANalytical B.V.
Lelyweg 1, 7602 EA Almelo
P.O. Box 13, 7600 AA Almelo
The Netherlands

T +31 546 534 444
F +31 546 534 598

info@panalytical.com
www.panalytical.com

Regional sales offices Americas
T +1 508 647 1100
F +1 508 647 1115

Europe, Middle East, Africa
T +31 546 834 444
F +31 546 834 969

Asia Pacific
T +65 6741 2868
F +65 6741 2166