



## Quality Control of $\text{Li}_2\text{CO}_3$ from Salars by XRD analysis using ARL X'TRA Companion benchtop XRD

### Authors

Dr. Simon Welzmler,  
Application Specialist XRD,  
Raphael Yerly,  
Product Manager XRD



Figure 1. ARL X'TRA Companion diffraction system

### Introduction

This application note presents the X-ray diffraction (XRD) analysis of lithium carbonate ( $\text{Li}_2\text{CO}_3$ ) from salars. Salars are salt flats or salt pans where lithium is extracted as  $\text{Li}_2\text{CO}_3$  or  $\text{LiOH}$  from brines or saltwater through a series of evaporation steps. Lithium carbonate is widely used in the production of batteries, ceramics, and glass, among others. Its purity and crystallinity are crucial to ensure the quality of the final product.

XRD is a non-destructive technique that provides structural information about crystalline materials. It is commonly used to identify and quantify the phases present in a sample and to determine their crystalline structure.

The results of XRD analysis provide valuable information about the quality and purity of  $\text{Li}_2\text{CO}_3$  from salars, which is essential for its application in various industries.

### Instrument

The Thermo Scientific™ ARL™ X'TRA Companion (c.f. Figure 1) is a simple, easy-to-use bench top XRD system for process control and more advanced applications.

The ARL X'TRA Companion uses a  $\theta/\theta$  goniometer (160 mm radius) in Bragg-Brentano geometry and a 600 W X-ray source (Cu or Co). The radial and axial collimation of the beam is controlled by divergence and Soller slits, while air scattering is reduced by a variable beam knife. An integrated water chiller is available as an option.

Due to the state-of-the-art solid state pixel detector (55x55 $\mu\text{m}$  pitch) the ARL X'TRA Companion provides very fast data collection and comes with one-click Rietveld quantification capabilities and automated result transmission to a LIMS.

### Experimental

Two powdered samples of  $\text{Li}_2\text{CO}_3$  from salars were measured in reflection geometry using an ARL X'TRA Companion with Cu  $K\alpha$  radiation (Ni filter). Each sample was measured for 10 minutes. Phase quantification was performed with Profex<sup>1</sup> (BGMN algorithm) using a fundamental parameters approach.

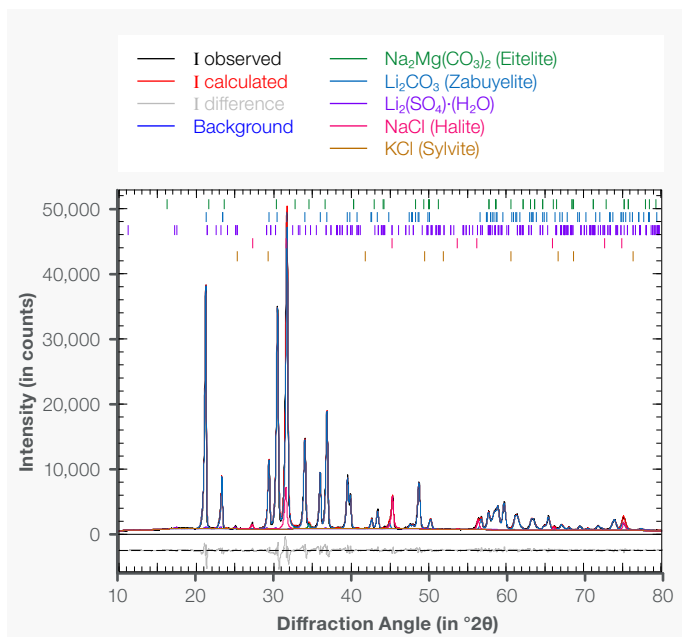


Figure 2. Rietveld fit of  $\text{Li}_2\text{CO}_3$  sample A (10 minutes)

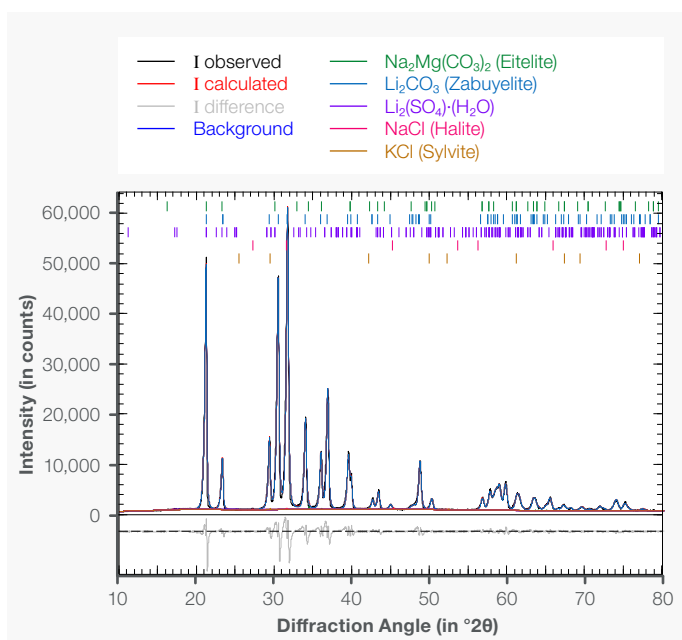


Figure 3. Rietveld fit of  $\text{Li}_2\text{CO}_3$  sample B (10 minutes)

Phase	Sample A (in wt%)	Sample B (in wt%)
$\text{Na}_2\text{Mg}(\text{CO}_3)_2$ (Eitelite)	0.7 (1)	0.2 (1)
$\text{Li}_2\text{CO}_3$ (Zabuyelite)	95.3 (2)	99.5 (2)
$\text{Li}_2(\text{SO}_4)\cdot\text{H}_2\text{O}$	0.7(1)	0.2 (1)
NaCl (Halite)	3.3(1)	0.1 (1)
KCl (Sylvite)	0	0
<b>Total Li (XRD/Reference)</b>	<b>18.0 / 17.2</b>	<b>18.7 / 18.5</b>

Table 1. Results of Rietveld refinement on  $\text{Li}_2\text{CO}_3$  samples (3 $\sigma$  in brackets)

## Results

In Figure 2 and 3 the refinements of two  $\text{Li}_2\text{CO}_3$  samples are shown.

Table 1 shows results from the refinements. The two samples are of different quality. Ascertainment of quality is important because phase-pure material with high total Li content is essential for further processing. The Li content can be indirectly ascertained from Rietveld refinements assuming stoichiometric phases with known Li content. Sample A is contaminated with 3.3 wt% NaCl and other minor phases, which results in a total Li amount of 18.0 wt%; this is in good agreement with the reference value of 17.2 wt% from chemical analysis. Sample B is of better quality and contains minor contaminations, which results in a total Li amount of 18.7 wt% (reference: 18.5 wt%). Both samples are completely crystalline.

## Conclusion

XRD is a powerful technique to assess the quality and purity of  $\text{Li}_2\text{CO}_3$  because it can indirectly measure the total Li content of a sample as well as contaminants which could impact the price and further processing of the material. The ARL X'TRA Companion XRD can collect data on  $\text{Li}_2\text{CO}_3$  in only 10 minutes, which allows the identification of contaminants and the quantification of the whole phase content. The one-click Rietveld refinement based on a fundamental parameter approach is an extremely robust method with high reproducibility. Thanks to these qualities, the ARL X'TRA Companion is the perfect solution for any process control or quality control task.

## Reference

1. N. Döbelin, R. Kleeberg, *J. Appl. Crystallogr.* **2015**, *48*, 1573-1580.

Learn more at [thermofisher.com/xtra](https://thermofisher.com/xtra)

**thermo**scientific

For research use only. Not for use in diagnostic procedures. For current certifications, visit [thermofisher.com/certifications](https://thermofisher.com/certifications)

© 2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. AN41502-EN-05-2023