

## Fast Ultrapure Water System with Inventive Bagtank Simplifies Process in Battery Excellence Lab

### Customer Challenge

The customer needed versatile balances and lab water systems to set up a new Battery Excellence Lab. They had the following criteria:

- Open connection to a new LIMS (not purchased yet)
- Balances for a broad range of laboratory analytics
- Fast access to ultrapure water for ICP-OES and ion chromatography applications

### Solution Provided

The customer attended an on-demand webinar and asked to be contacted by an expert. Together with a Sartorius Virtual Account Manager, they discussed the customer's sample preparation workflow and criteria for semi-micro balances and ultrapure water.

Balances are involved in several steps of the battery workflow. These include pre-analytical preparation steps and material characterization, such as calibration curves preparation, density determination, specific surface area, Karl Fischer, and electrode grammage.

Sartorius recommended the Cubis® II semi-microbalance (capacity 220 g, scale interval 10 µg) with an advanced user interface MCA, automated draft shield with built-in ionizer, Qapp utilities package QP3, and foot switch door opener. The customer liked these features and was especially interested in the data integrity, electronic signature, user password, and open LDAP interface to connect with a LIMS they plan to acquire.

The customer also required 2 dispensing of 5 L of ICP- and chromatography-grade lab water per day. They requested a lab water system that delivers prompt access to water after regeneration or filling of internal tanks and is easy to maintain. Sartorius recommended the Arium Pro UV 18,2 MΩ.cm-1 with its unique Bagtank, storage of pre-treated water. The disposable bags allow fast and easy replacement with a new bag instead of the time-consuming cleanings for tanked systems.

### At a Glance

- Future-proof, open interface balance with data security
- Qapp package includes grammage determination YAPP305
- Easy to maintain lab water system with instant access to ultrapure water

For more information, visit:



### Case Profile

**Company Type:**  
European Company

**Industry:**  
Battery Excellence Lab

**Company Size:**  
< 50 employees

### Outcome

The customer selected both Cubis® II and Arium® Pro UV for their sample preparation and characterization workflows. The local Sartorius Service Team installed the balance and the lab water system, and conducted the necessary user training for cartridge replacement and other routine maintenance. Discussions are underway about a yearly supply of ash-free filter paper.