

Built-In Ionizer Enables Reliable Weighing in Battery Material Extraction Workflow

Customer Challenge

Customer specializing in battery material extraction needed balances to equip their R&D and QC labs. They had the following criteria:

- Hassle-free weighing of static powders from lithium salts extraction
- Accurate and reliable weighing before chemical and physical characterization
- Easy collection of weighing data for further Design of Experiments analytics (DoE)

Solution Provided

The customer first contacted a Sartorius Salesteam member to ask about possible solutions to improve their weighing workflow.

After completing the Lab Weighing Academy training, they learned about the factors that impact weighing accuracy and identified statics and poor leveling as their primary challenges. Next, the customer requested on-site testing of different balance models.

Sartorius arranged for an on-site demonstration of the Cubis® II balance. This balance offers five-digit precision, an advanced user interface MCA, built-in ionizer, motorized leveling and QP2 onboarded Qapp software. During the trial, the customer found that the stability, reproducibility, easy cleaning and data transfer met all their criteria. Most importantly, the motorized auto levelling and built-in ionizer addressed electrostatic disturbances.

For less demanding applications, like Check Weighing, the customer liked the Entris® II Essential, with two-digit precision, internal calibration isoCAL, a large 182x182 mm weighing pan, fast stabilization time, and high chemical resistance.

At a Glance

- Built-in ionizing technology reduces electrostatic effects
- Automatic and motorized leveling
- Open interface for accessible transfer of data

Cubis® II Videos:

For more information, visit:



Case Profile

Industry:
Material Extraction

Company Size:
100 000+ t/year Lithium salts

Outcome

The customer selected both Cubis® II and Entris® II models for their R&D, Quality Control Labs, and off-line production for non-destructive controls. The local Sartorius Service Team installed and calibrated the balances, and provided training on routine cleaning and maintenance. To secure balance longevity and reduce downtime, a yearly maintenance contract was signed.