

Next-Gen Battery Innovation Lab Invests in Fully-Automated Weighing Technology

The customer is an innovation lab in the battery space and is heavily invested in developing tomorrow's solid-state batteries, as a safer, lighter and more compact alternative to current offerings.

Testing Challenge

- Necessity for precise weighing within a dry and inert atmosphere to prevent external variables from influencing results
- Requirement for hands-free operation to maintain workflow efficiency and minimize contamination risks
- Need for an effortless method to capture and analyze weighing data for further Design of Experiments analytics (DoE)
- Fast and precise way to measure moisture content in plastic tubes components of the Battery Management System (BMS)



The Solution

- Cubis® II semi-micro balance featuring an automatic draft shield, a built-in ionizer to neutralize static, and motorized leveling for optimal accuracy. The balance was combined with a footswitch command
- Each weighing datapoint was openly collected via RJ45 for further Data Analytics
- MA160 Moisture Analyzer (capacity 200 g; moisture determination > 1%) backed by its Method Development Assistant function



For more information,
visit www.sartorius.com

The Value

- Accurate, traceable and reliable weighing data even in a challenging environment
- Local Sartorius Service Team provided comprehensive support, including installation, calibration, training, and a yearly maintenance contract, ensuring sustainable longevity of the balances and minimizing potential downtime
- As molding temperature depends on the moisture content of in-coming goods, fast moisture determination helped for instant machine parameter settings

SARTORIUS