

Cubis® II High-Capacity Micro Balances for On-Body Self-Medication

The Company's on-body injector simplifies treatment, reduces healthcare costs, enables the administration of new therapeutics, and improves patient experience. It facilitates versatile subcutaneous delivery of up to several milliliters for three-day wear and offers programmable, connected medication administration outside of hospital settings.

Testing Challenge

- Stable measurement of the bolus
- Very brief shot intervals (seconds) over a long period of time (days)
- Electrostatic charges due to plastic components

The Solution

- Outstanding Cubis® II High-Capacity 61 g microbalance that ensures reliable flow and accuracy (F&A) determination over several days
- Integrated deionizer paired with coated draft-shield

The Value

- Optimized shot pattern during uninterrupted testing periods
- Reduced need for frequent weighing reservoir emptying
- Electrostatic charge removal for a stable acquisition of weigh value



For more information,
visit www.sartorius.com

SARTORIUS