

Cubis® MPS Balances for Insulin Pump Testing

The Company specializes in manufacturing wearable insulin patches. Its wearable pump is designed to be fully disposable and has a capacity to hold a sufficient amount of insulin for up to 3.5 days. Engineered to administer both basal and bolus insulin doses as needed, the pump is also intended to connect wirelessly to a user's smartphone, thereby simplifying the management of insulin dosages.

Testing Challenge

- Basal micro-flow rate and bolus dose accuracy determination
- Over a period of up to 7 days
- The calibration lab, situated adjacent to production, is susceptible to vibrations



The Solution

- Cubis® MPS 0.01 mg balance designed for volume calibration, with large reservoir, evaporation trap, and automated draft shield opening
- Weighing system is isolated from the balance interface and the entire test setup is protected from external disturbances by a protective box

The Value

- Hands-free collection of experiment data
- Minimal to no impact from environmental fluctuations
- Dynamic conversion of weighing value into long-term flow-rate evaluation



For more information,
visit www.sartorius.com

SARTORIUS