

Cubis® Balances for Infusion Systems Development

The Company is engaged in the development, production, and distribution of dialyzers and automated peritoneal dialysis systems designed to treat immune disorders, infectious diseases, and trauma-related conditions. The Company's Research and Development division was established with the objective of providing internal IT solutions to support the Company's global business operations.

Testing Challenge

- Every piece of Equipment Under Test (EUT) is required to function uninterrupted and flawlessly for the entirety of the testing phase, which spans several months
- Broad range of nominal flow rates
- Compliance with medical device regulations



The Solution

- Cubis® lab balances with a capacity from 220 g to 20 kg
- Accurate and linear dynamic flow measurement
- Open access to balance data for easy acquisition of analytical parameters through a customized software
- Qualified installation with on-site IQ/OQ/PQ

The Value

- Stable data output is driven by consistent weighing performance, even in challenging environments
- Easy transfer of analytical method to other reference labs in the Company, supported by robust documentation
- Minimal drift ensured by monolithic weighing system and electrostatic charge removal



For more information,
visit www.sartorius.com

SARTORIUS