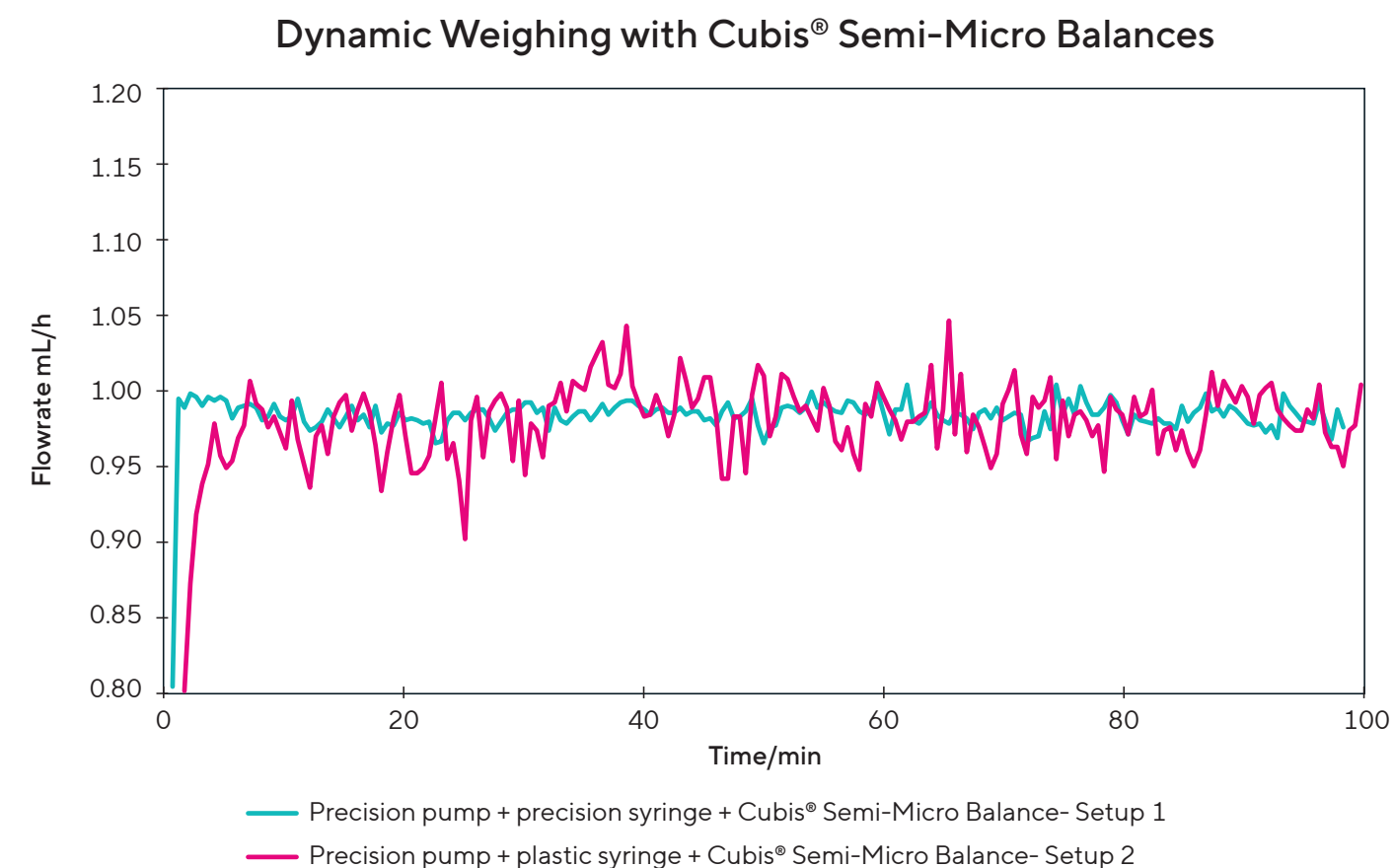


Cubis® Semi-Micro Balances to Assess Syringe Precision

The Company specializes in providing advanced automated infusion systems for patient care, ensuring high standards of safety and efficiency in medical therapies. The Company offers a comprehensive range of products and services, including clinical IT solutions and disposable products, to support healthcare professionals and enhance patient outcomes..

Testing Challenge

- Ensure that fluctuations in flow rate are not artifacts caused by balance drift
- Compare single use syringes with precision syringes to assess syringe accuracy and batch-to-batch homogeneity
- Demonstrate the accuracy of syringe-infusion pump combination over a certain period of time



The Solution

- Cubis® II Semi-Micro balance's for accuracy measurement in dynamic weighing
- Consistent flow rate near 1 mL/h, demonstrating compliance with ISO 7886-2 and IEC 60601-2-24 accuracy requirements
- Highly precise measurements, crucial for ensuring accurate medication dosing

The Value

- Disposable plastic syringe showed larger deviations, indicating the impact of accessories on measurement accuracy
- Reliable balance's performance in dynamic weighing
- Highlights the importance of using high-quality accessories to maintain accuracy



For more information,
visit www.sartorius.com

SARTORIUS