SARTURIUS

Application Guide

May, 2024

Keywords or phrases:

Cleanliness, particulate matter, moisture determination, residue of ignition, total solid content, density, specific gravity, filter weighing, biotherapeutics, chemicals, classical pharma, cosmetics, F&B, medical device, environmental labs, testing labs, automotive, battery manufacturer

More Sample Preparation

Fundamental PreAnalytics for Compliant Flexibility

Sartorius provides an extensive selection of lab essentials tailored for the initial stages of sample handling and preanalytical processes in quality assurance labs, catering to diverse characterization requirements. Our lineup includes accurate lab scales, high-purity water purification systems, sophisticated filtering equipment, and comfortable pipettes and tips. These instruments are engineered to support precise characterization of particulate matter, moisture levels, total solid content (TSC), residue of ignition, and density, even under challenging conditions. Additionally, they are proficient in preparing for detection of trace impurities like phthalates and PFAS in different substances, guaranteeing thorough and precise analytical outcomes.







Lab Balances

- Preinstalled Applications to Comply with Main Pharma Regulations
- Customizable and Upgradable Balances to Meet vour Actual and Future Needs

Pipettes and Tips

- Versatile Range of Mechanical and **Electronic Pipettes**
- Matching Tips for Accurate Results

Lab Water Systems

- Extensive Portfolio to Get the Right Lab Water Where you Need it
- Different Set-up to Adapt to your Requirements

Filtration & Retention

- Broad Choice of Membrane Types and Diameters
- Configurable Filtration Manifold. Support and Funnels

■ Lab Balances		Pipettes & Tips Lab Water Systems	Filtration Devices	
Analytics	Topic	Details	Sartorius Solution	Link
Inert Atmosphere, Glove Box	Optimizing Balance Performance in an Inert Gas Atmosphere	Application highlight. Precise weighing within an argon gas box, necessary conditions for maintaining accuracy and challenges posed by glove use and spatial constraints.	■ Cubis® II	母
Density 	Comprehensive Guide to Density Determination Using Lab Balances	Guide. Principles of density measurement, traditional and modern techniques, and role of lab balances in data accuracy for QC and R&D.	■ Entris® II ■ Cubis® II	⊕
Cleaning ■	•	Guide. Best practices for maintaining Cubis® II balances, daily and extensive cleaning procedures, chemical compatibility, and dedicated cleaning App.	■ Cubis® II	@
PFAS	PFAS: Environmental Persistence and Health Risks	Expert insight. Environmental presence and health risks of PFAS, research on their detection, effects, and passive sampling devices to assess their bioaccumulation.	■ Arium®	⊕
Electrostatic Charges	Stable weighing protected from outer static effects	d Flyer. Cubis® II balances featuring coated draft shields acting as Faraday cage, neutralizing static interference and ensuring precise measurements in lab settings.	■ Cubis® II	₩
Protein Concentration	Ultrafiltration in Clinical Diagnostics: A Pre- Validation Guide	Guide. Pre-validation process of GLU ultrafiltration devices in clinical sample prep, protein concentration and recovery for reliable diagnostics.	■ Vivapore® ■ Vivaspin®	⊕
Heavy Metals	Groundwater Sample Filtration for Heavy Metal Analysis	Application note. Sartolab® P20 Plus filter's efficient with groundwater samples for heavy metal testing, superior throughput and minimal metal ion release. In collaboration with Eurofins.	■ Sartolab® P20 Plus	₩

Germany

Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Strasse 20 37079 Goettingen Phone +49 551 308 0

For further contacts, visit www.sartorius.com

USA

Sartorius Corporation 565 Johnson Avenue Bohemia, NY 11716 Phone +1 631 254 4249 Toll-free +1 800 635 2906