



Spotlight on Our Lab Essentials


2025 Edition

Simplifying Progress

SARTORIUS

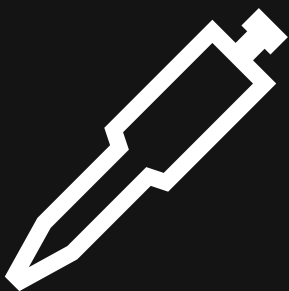
The Latest Solutions for Your Lab

Working in a scientific lab environment means you face growing challenges every day. At Sartorius, our daily mission is to help you enhance lab efficiency. That’s why we continuously develop innovative products and services to facilitate and accelerate your workflows. Find out how you will benefit from our latest lab solutions and our longstanding expertise.

 **Find out more**
For more information, please visit
www.sartorius.com



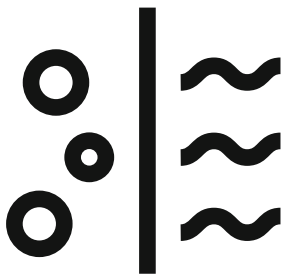
Lab Weighing



Pipetting & Dispensing



Lab Water



Lab Filtration
Reagents & Consumables



Microbiological Analysis



Software

Did you know?

Sartorius is committed to the long-term goal of fully decarbonizing its business activities by 2045. The company plans to achieve climate neutrality by continuously decarbonizing along the value chain and offsetting unavoidable residual emissions in line with the Paris Climate Agreement.

We aim to fully decarbonize our business activities by 2045.

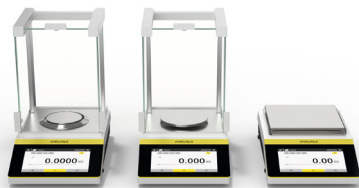
2030 mid-term targets

~10%	Zero	100%
average reduction of CO ₂ eq emission intensity p.a. (Scopes 1-3)	avoidable emissions in Scopes 1 and 2	electricity from renewable sources
<hr/>		
>75%	Zero	
revenue with products designed according to circularity principles	operational waste to landfill	

2045 long-term target

Net-zero emissions

Decarbonization of the entire supply chain in collaboration with suppliers and customers



Quintix® Pro Balances are manufactured using 100% renewable energy. Equipped with energy efficient designs such as automatic dim light to conserve power when not in use.

Picus® 2 Electronic Pipettes and Pipette Tips are manufactured using 100% renewable energy. Production waste is less than 5% of total plastics used.



Sartorius Refill Tower earned ACT label in 2024.

Download Sustainability Fact Sheets for our products here!

Sustainability fact sheets are concise documents highlighting key information about an organization’s sustainability practices, initiatives, and impacts. They include data and statistics on environmental, social, and economic efforts, serving as a tool for transparency and accountability to stakeholders.





Entris® II Basic Balance

Best Value in Its Class for Basic Weighing Tasks

No matter what you're weighing, our Entris® II balance is always the right choice. Our Entris® II Essential line is the only balance in its class offering isoCAL, LED touch technology, and 12 built-in applications at a budget price. With over 40 models from which to choose, you will surely find a model to meet your basic weighing needs.

Features and Benefits

- High chemical resistance – ensured using parts made from hard-wearing Polybutylene terephthalate (PBT), stainless steel and glass
- isoCAL – automatic internal calibration and adjustment
- Fastest stabilization time in its class
- Real „PC-direct“ feature
- 12 built-in applications with GLP | GMP compliant printout | data output

Entris® II Advanced Balance

Every Sartorius balance offers quality, value, and consistency

Our Entris® II Advanced line gives you additional value to the Entris® II portfolio, with over 38 models offering benefits such as real-time level support, integrated protection systems, CalAuditTrail, a graphic touch display and 13 built-in application programs.

Features and Benefits

- Simplifies balance leveling with built-in real-time level support
- Increases reliability of weighing results using integrated protection systems
- Controlled access to balance settings with user management
- Gapless documentation thanks to CalAuditTrail
- Even easier usability with graphic touch display
- Provides gapless documentation with automatic documentation of calibration and levelling processes, and date and time stamp
- One additional built-in application: Pipette Smart Test



Cubis® II Premium Balance

Data Integrity and Connectivity

Need to comply with 21CFR11? Need the fastest stability on the market? Cubis® II – Your Balance, Your Way.

Features and Benefits

- The world's highest resolution in a micro weighing instrument with up to 60 million divisions
- Integrated Software - no need for additional computer
- Full end-to-end data integrity. 21 CFR Part 11 compliance, integrated audit-trail, state-of-the-art user management
- Error-free operation. Individual QApp workflows, motorized auto-leveling*
- Integrated climate sensors
- Individual sample holders
- Outstanding service support. Integrated status center, integrated service functions preventive maintenance on accredited standards
- **NEW:** Cubis® II Ultra High-Resolution High-Capacity and Semi-Micro balances which guarantee the highest level of precision with each measurement in everyday conditions.

* For all models up to a maximum capacity of 8.2 kg

Quintix® Pro

Redefining the Standard

It's easy to overlook the standard lab balance. To expect the ordinary. Until something new comes along that redefines what a standard balance can be.

Quintix® Pro Laboratory Balances deliver the performance you need, with added premium features that raise the bar in usability, flexibility, and eco-conscience design.

Features and Benefits

- Enhance the user experience with new premium features
- Offering a wide range of use in different environments and applications
- Energy efficiency and reduced carbon footprint minimize environmental impact from the manufacturing plant to the bench



Efficient
Product Use



Climate-
conscience
Production



Sustainable
Packaging





Tacta® Mechanical Pipette

Features & Benefits

- Comfortable to hold due to the ergonomically designed handle
- Low pipetting and tip ejection forces that reduce the risk of Work Related Upper Limb Disorder (WRULD)
- Controlled and smooth tip ejection with the new Sartorius Optiject technology
- Calibration adjustment to provide accurate results for various liquid types
- Easy to clean, with only three parts to disassemble



Sartorius Pipetting App

A must-have for all Picus® 2 pipette users, the Sartorius Pipetting app is your handy companion for sample preparation workflows and pipette management. To stay up to date, use the Pipetting app to update your Picus® 2 pipettes with the latest features. Whether you are a seasoned scientist or a novice in the field, this app offers accuracy and reproducibility for your liquid handling workflows.

The app features an ever-growing database of configurable protocols that can be adapted according to your specific application. The app provides step-by-step guidance for various workflows, and when used with Picus® 2 pipettes, it automatically selects the correct settings. Complete your workflow using Picus® 2 pipettes, other electronic pipettes, or any mechanical pipette and export the logfile for your records.

Visit www.sartorius.com for updated list of workflows.

Volume Adjustment

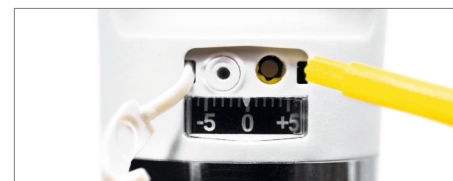
Use the Sartorius Optilock feature to prevent accidental volume changes during pipetting. Either hold down the volume lock button with one hand and adjust the volume with the other, then release to lock, or with just one hand, slide the volume lock up, adjust the volume and slide down to lock again. The choice is in your hands.

Large, Clear Display

The large, easy-to-read display helps you to see all four digits of the set volume from a distance without straining your eyes. The volume is easy to read even when the pipette is angled – eliminating the need to turn your head into an uncomfortable position.

Calibration Adjustment

Tacta® is very easy to adjust, for various types of liquids, using a simple adjustment key. The adjustment scale shows the degree of adjustment. By noting this value for a specific liquid, you can return to that setting any time. Easy calibration adjustment ensures accurate results with a variety of liquids.



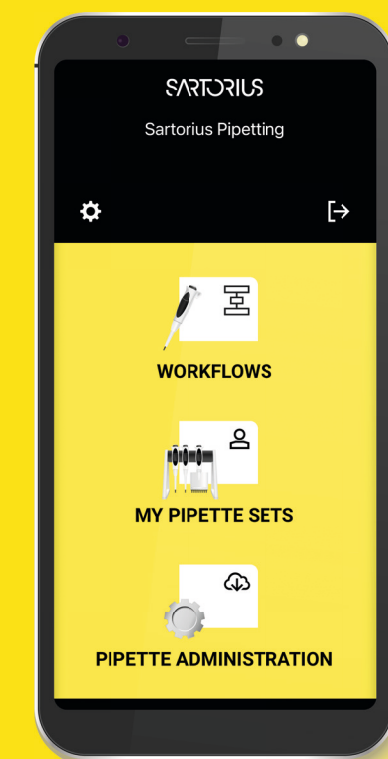
New Picus® 2 Electronic Pipette

The Most Sophisticated and Ergonomic Pipette Ever!

The Picus® 2 pipettes ensure reliable and repeatable pipetting results with unbeatable ergonomic design that is kind to your hand. Picus® 2 is the tool of choice for everybody from a student to experienced laboratory professional. It is as intuitive to use as a mechanical pipette and yet offers a broad range of options even for the most advanced users.

Features and Benefits

- Highest level of ergonomics provided by the uniquely low weight, light electronic tip ejection and comfortable handle design
- Extensive range of pipetting modes reduces the needed pipetting steps and speeds up work
- Electronic brake and piston control system provide outstanding accuracy and repeatability of pipetting results, independent of the user
- Intuitive user interface with language options, enables ease of use
- Adjustment wheel offers extremely fast volume setting and menu navigation
- Optiload enables perfect tip sealing for accurate delivery from each channel



Download the App:



Sartorius Optifit® Tips and Safetyspace® Filter Tips are designed to fit perfectly Sartorius pipettes for maximum pipetting accuracy and reproducibility.

Features

- Perfect and soft sealing with Sartorius pipettes
- Unique safety space between filter and sample
- Extremely even and durable surface in low retention tips
- Packing options for every lab and color-coding to match your pipette

Benefits

- Ensures precise and repeatable results of your pipette
- Protect sample purity with Safetyspace® filter tips and purity certified Optifit® tips



Sustainable Packaging



Optimized Transport & Logistics



Optimized Production

Arium® Mini Family

Maximum Flexibility in One Compact Design: Lab Water Purification with Arium® Mini and Arium® Mini Extend

The compact Arium® Mini water purification systems have been designed to meet your Type 1 ultrapure water needs, handling up to 10 liters per day. Developed according to Type 1 water requirements for general, life-science and analytical applications, this system guarantees reproducible results and streamlined processes. Operate effortlessly using the large, touch-activated screen and user-friendly menus. Choose from the Arium® Mini or Arium® Mini Essential if using pretreated water feed, or opt for the Arium® Mini Plus with innovative bagtank technology for a simple connection to your potable tap water supply.

Introducing the newest member of the Mini family: Arium® Mini Extend. Experience maximum convenience with the flexible, removable dispensing unit, keeping you ahead in your lab routines. Explore our extensive range of Arium® Mini lab water systems and choose the one that best fits your needs.



Features and Benefits

- Flexible: Different versions for different inlet water and connection options, as well as the desired water withdrawal
- Compact: Space-saving with a width of only 28 cm (+ 9 cm for versions with attached dispensing unit)
- Intuitive: Color touch display, with direct access to all important information and dispensing options
- Innovative: Depending on the type of system, our unique Bagtank technology saves time as no intensive tank cleaning with hazardous substances is required
- Reliable: Delivers consistently high water quality for reliable and reproducible results

Arium® Pro

Tailored to Your Specific Requirements

As a reliable source of ultrapure water, the Arium® Pro series offers a flexible and modular system which, compared to conventional devices, demonstrates excellent added value. All systems meet and exceed the ASTM Type 1 water quality standards and ensure the best reproducible results in their class. Up to 2 liters of consistently high-quality ultrapure water with a conductivity of 0.055 µS/cm can be dispensed per minute. The ultrapure water is virtually microorganism-free when an Arium® SterilePlus final filter is used. The patented Sartorius technology, the SD card slot, the long service life and low maintenance requirement distinguish the Arium® Pro systems as an easy-to-use, economical and reliable Type 1 ultrapure water system.

Features and Benefits

- Modular – System selection specifically for your application
- Flexible - Perfect integration into any laboratory
- Easy to use – Display with touch function and intuitive menu
- Fast – Favorites function with direct access for recurring volumes

Choose Your Smart Station

Connect up to 3 Smart Stations to your ultrapure water system for maximum flexibility. Supply ultrapure water from Arium® Pro and Arium® Comfort.



Arium® Comfort

Space-saving Twin Technology

Arium® Comfort I is an easy to use system for producing ASTM Type 1 ultrapure water and Type 3 pure water. The system contains state-of-the-art reverse osmosis technology and a unique cartridge specifically for the production of the highest ultrapure water quality.

Compared to conventional water systems, the Arium® Comfort I optimizes water consumption using the integrated iJust control unit. This unique touch display with intuitive menu navigation ensures the utmost ease of use.

With the optionally integrated TOC monitor, its compact design, the flexible display and the SD card slot, the Arium® Comfort I is the ideal choice for demanding laboratory applications.

Features and Benefits

- Time-saving – Use of innovative bag technology, eliminates costly tank cleaning
- Optimized water consumption – automatic with iJust
- Easy to use – Display with touch function - and intuitive menu
- Quick – Favorites function with direct access for recurring volumes

Feed Water	Arium® Mini	Arium® Mini Plus	Arium® Mini Essential	Arium® Mini Plus Extend	Arium® Mini Essential Extend
Tap Water		✓		✓	
Pretreated water (RO DI line)			✓		✓
Pretreated water (from supply container)	✓				
Flexible Removable Dispenser			✓		✓
Bagtank Technology	✓	✓		✓	
Purified Water					
Type 1	✓	✓	✓	✓	✓
Type 3		✓		✓	



Sartolab® RF | BT Vacuum Filtration Units

Sartolab® RF | BT vacuum filtration units are convenient filtration units designed for the filtration of volumes from 50 mL to 1 L. Sartolab® RF as a complete system includes a receiver flask or a conical tube to the filtration funnel. Sartolab® BT is a bottle top filter (filtration funnel) without a receiver flask, enabling customers to use their own receiver flasks and | or to expand the filtration capacity, depending on the particle load of the filtered liquid, by filling more than one receiver flask.

All filtration units can be used as stand-alone system or with the Sartolab® Multistation, which allows parallel filtration of up to 6 samples.



Sustainable Packaging



Optimized Transport & Logistics



Optimized Production

Sartoclear® Dynamics Lab

Quick and Efficient Harvest of Animal Cell Cultures

Sartoclear® Dynamics Lab ready-to-use-kits combine a filter aid with a vacuum filtration unit for clarification and sterile filtration in a single step of high density mammalian cell cultures.

Features

- Cell culture filtration of up to 1L in single step
- Filtration unit with 0.22 µm PES membrane
- Vacuum filtration unit with filter aid included in single kit
- Sartoclear Dynamic Lab kits compatible with Sartolab® Multistation

Benefits

- Fast result with high flow rates and single step filtration
- Efficient - No need of centrifugation step
- Hands-free parallel filtration with the Sartolab® Multistation without need of extra vacuum connector installation and time-consuming stabilization

Minisart® Syringe Filters

No. 1 Choice for Your Filtration Applications

Innovative design coupled with the largest surface areas and fastest flow rates make Minisart® syringe filters the ideal choice for all your syringe filtration needs.

Features

- Large effective filtration area (EFD) of 6.2 cm²
- Available in wide range of configurations
- Low adsorption
- Superior flow rate

Benefits

- Fast simple, and economical
- Low hold-up volumes and maximal recoveries
- Low adsorption of analytes
- Maximum chemical compatibility
- Minimum extractables or leachables
- Bidirectional use possible
- 100% optical integrity test
- Membranes with a wide choice of pore sizes
- HPLC Certification available

Alternative Tip

Are you looking for a high-throughput, easy to use syringeless system for your HPLC sample prep? Check out our Claristep® Filtration System!



Sartolab® Multistation for Parallel Filtration of up to 6 Samples

Sartolab® Multistation is a stand uniquely designed to hold one to six vacuum filtration units, allowing simultaneous filtration of up to six samples. Easily install your vacuum filtration units in the Multistation for quick and easy filtration without the need for extra connectors and time-consuming stabilization.

Sartolab® Multistation works with Sartolab® RF|BT vacuum filtration units for filtration of 50 mL to 1 L samples.



Sartolab® P20 Pressure Filtration Devices

Sartolab® P20 filtration devices are ready-to-use filters for the clarification and sterile filtration of media and aqueous solutions in batches from 100 mL to 10 L. They are available either with 0.2 µm or 0.45 µm PES membranes, with or without a prefilter made of high purity quartz microfibers. An additional version containing the quartz microfiber prefilter only is also available for clarification purposes.



Quantitative (Ash-Free) Filter Papers

In quantitative analyses, either the filter cake or the filtrate is analyzed.

In gravimetric analyses, the filter cake, along with the filter paper, is calcined to determine the ash content. Filter papers with an ash content of less than 0.01% are considered ash-free and can be used for such analyses.

Ash-free filter papers are also essential when analyzing the filtrate to avoid contamination from foreign substances. These papers are exceptionally pure, made from cotton linters, and acid-washed during production to remove trace elements.

- Applications**
- Gravimetric and quantitative analyses

- Properties**
- Made of cotton linters with over 98 % alpha-cellulose content
 - Ash content < 0.01 %
 - Wet strength
 - 6 Grades for particle retention from approx. 1 µm to 15 µm



Qualitative Filter Papers

Qualitative filter papers are used primarily for general laboratory applications where the primary goal is to separate solids from liquids without the need for precise measurement of the retained material. The filtrate only will be analysed.

- Applications**
- General laboratory tasks (separating solids from liquids, clarifying, removing particle matter)

- Properties**
- Plain or creped filters
 - Ash content < 0.1
 - Made of cellulose or refined pulp and cotton linters with over 95 % alpha-cellulose content
 - Wet strengthened



Glass & Quartz Microfilter Filters

They are essential for applications requiring higher temperature resistance, fine particle retention, and chemical resistance. They are available either with binder or free from any additives.

- Made from borosilicate glass fibers:**
- Glass microfiber filter with binders
 - Glass microfiber filter without binders

- Made from high-purity quartz fibers:**
- Quartz microfiber filters
 - Tempered microfiber filters

- Properties**
- High flow rate for fine particles
 - Higher chemical and temperature resistance compared to filter papers
 - Mechanically more stable than filter papers



Extraction Thimbles

Our extraction thimbles have been designed to be compatible with the Soxhlet apparatus, making them ideal for a variety of applications. These include environmental monitoring tasks such as the separation of particulates like dust and aerosols, as well as the analysis of gas or air streams.

- Composed of high-purity cellulose and glass fiber for uncontaminated results
- Engineered for superior mechanical stability to withstand rigorous use
- Glass fiber thimbles offer remarkable resistance to high temperatures
- Designed for a precise fit across a diverse range of extractors
- Exhibits outstanding chemical resistance, ensuring reliability in various extraction processes.



Sterisart® Universal | Gen 4

Where Convenience Meets Compliance

The Sterisart® Gen 4 Pump features a compact, universal design that easily fits into various workspaces, including benchtops in laminar airflow cabinets, RABS, and sterility testing isolators. Select the model that meets your needs with the convenience of built-in upgradability.

Features and Benefits

- Universal sterility testing pump designed for benchtop use or for integration in isolators
- Low profile and compact footprint
- Flexible, modular and open design
- 316L mirror finish stainless steel
- Sleek design for impeccable hygiene: easy disassembly and smooth contours
- Real-Time pressure detection in the Sterility testing canisters
- 7" high performance touch screen display with integrated barcode scanner
- Built for 21 CFR Part 11 & EU GMP Annex 11: Advanced software for sterility testing featuring user access control, audit trail, secure data handling, electronic signatures, and reports

Sterisart® NF System

Your Choice For Sterility Testing

The Sterisart® NF System is a unique and completely disposable technology based on the membrane filtration method used in conjunction with the Sterisart® Universal Pump which allows for the sample and media transfer within the Sterisart® NF System. Both work together to enable the sterility testing of pharmaceutical & medical device products in full compliance with international pharmacopeias.

Features and Benefits

- Unique Septum for Aseptic Sampling
- Pre-installed color-coded tube clamps for easy and reliable assignment to each Sterisart® NF container
- Product/lot number identification for traceability of the product for documentation and security
- Modular design pump which is easy to decontaminate

Microsart @filter & Microsart @media

Advancing Membrane Transfer for Microbial Enumeration

Microsart® @filter combined with Microsart® @media offer a unique hands free membrane transfer technique minimizing the risk of contamination during your microbial enumeration test.

Features and Benefits

- The innovative membrane transfer is a touchless transfer to the agar avoiding the introduction of air bubbles ensuring perfect membrane to agar contact
- The design of the Microsart® @filter base fits perfectly onto the Microsart® @media for reliable transfer of the membrane
- The reliable membrane transfer reduces handling and minimizes sources of contamination resulting in more reliable testing results

MD8 Airport – Portable, Reliable, Fast

MD8 Airport is the air sampler for the pharmaceutical industry, biotechnology, food and beverage industry, for hospitals, and for measurements within the scope of environmental protection and occupational safety.

Features and Benefits

- Battery operated and portable – so it can be used everywhere
- Battery power level clearly indicated so constant performance during sampling is guaranteed
- Ergonomically designed and easy to clean
- User friendly prompting with the option of four languages: English, French, German and Spanish
- Flexible options for customized use including adjustable air throughput and air flow rate
- Parameters last used are stored even after automatic shut-off
- Can be calibrated on-site
- Utilizes Gelatin Membrane Filter Method, allowing for reliable and accurate results





Sartobind® Lab IEX

Fast, Flexible and Future-Proof Charge-Based Separation

Sartobind® IEX Lab incorporates the state-of-the-art in ion exchange chromatography (IEX) membranes, enabling flexible biomolecule purification that is both fast and scalable. It is ideal for protein, virus and nucleic acid research, as well as the development of downstream processes to purify novel therapies and vaccines before scaling up to clinical and commercial production.

Features and Benefits

- **The Fastest Matrix:** Full speed ahead. Rapid flow rates and reduced residence times take you from equilibration to pure molecule in a matter of seconds, not minutes.
- **Productivity Boost:** Push the limits. Macroporous membranes with high ligand densities deliver maximum yields of even the largest viruses without size exclusion effects.
- **Plug and Play:** Collapse your workflow. Ready-for-use units avoid the need for column packing and buffer degassing, while enabling convenient inline prefiltration.
- **Process-Ready Technology:** Seamless scalability. Future-proof by design, ensuring ease-of-transfer for your most promising targets to clinical and commercial production.

Sartobind® Rapid A Lab

To Affinity and Beyond

Sartobind® Rapid A Lab incorporates the state-of-the-art in membrane affinity chromatography (AC) technology, offering fast, flexible and future-proof antibody purification. It is an ideal solution for antibody research, as well as downstream process development for novel biologics before starting clinical and commercial production

Features and Benefits

- **The Fastest Matrix:** Full speed ahead. High flow rates and reduced residence times take you from equilibration to pure antibody in less than two minutes.
- **Productivity Boost:** Push the limits. Maximum dynamic binding capacity supports the preparation of your antibody sample to the highest possible yield and purity.
- **Plug and Play:** Collapse your workflow. Ready-for-use units avoid the need for column packing and buffer degassing, while enabling convenient in-line prefiltration.
- **Flexibility Built In:** Accept no limitations. Enhanced pressure resistance in a platform that ensures reliable purification with or without a liquid chromatography system.
- **Process-Ready Technology:** Seamless scalability. Future-proof by design, ensuring ease-of-transfer for your most promising targets to clinical and commercial production.

Vivaspin® Centrifugal Concentrators

Flexible and Efficient Protein Concentration

Vivaspin® products offer a comprehensive range of devices for fast protein concentration from 0.5 ml to 20 ml volume capacities. The large range of different molecular weight cut offs, from 2 kDa to 1000 kDa and 0.2µm offer a solution to every concentration need. Unique Vivaspin® 20 can be pressurized and centrifuged simultaneously for processing difficult samples.

Features and Benefits

- **Guaranteed sample recovery** – no risk of concentrating to dryness
- **Vertical membrane configuration**
- **Easy to find the perfect ultrafiltration device** for your application due to large product selection. Widest volume capacity range and molecular weight cut off ranges
- **Important:** Centrisart 1 and Vivacell 100 have recently joined the Vivaspin family as Vivaspin® Filtrate and Vivaspin® 100, respectively

Vivaflow® SU Crossflow Devices

Intuitive, Effortless and Sustainable TFF for the Lab

Vivaflow® SU is a next generation plug-and-play tangential flow filtration (TFF) cassette, for ultrafiltration and diafiltration of intermediate sample volumes. It has been specifically designed for ease of use in research and development laboratories.

Features and Benefits

- **True Plug and Play** - Link the cassettes quickly and securely to a suitable pump and reservoirs, right out of the box
- **No Added Costs** - All-in-one design avoids the need for expensive cassette holders and TFF systems
- **Performance Optimized** - Advanced flow path and fixed occlusion restrictor eliminate the effort that is usually needed for TFF
- **Maintain Sample Integrity** - Improve safety for you and your sample with no need for flushing or cleaning



Sustainable Packaging



Optimized Production



Ingenix Suite - Ingenuity in the Lab



Simplifying Your Lab Workflows: Elevate Accuracy, Efficiency, and Compliance for Pipettes and Lab Balances Seamlessly

Ingenix Suite eliminates manual steps by centralizing and automating data transfer, job management and reports generated by your instruments (balances, pipettes). Increase your lab efficiency, data quality and enhance compliance.

With its flexibility and openness, Ingenix Suite works with or without an existing ELN/LIMS system on site. Because your lab needs will never be covered by a single vendor, Ingenix Suite is open (module dependent).

Ingenix Suite Balance Device Manager Module

- Manage your entire Cubis®II MCA balance device across all labs with unlimited connections
- Centralize weighing data in an SQL database for easy access
- Easily update Cubis® II MCA firmware remotely
- Enhance service by sharing diagnostic information remotely

Ingenix Suite Pipette Check Module

- Quickly verify pipette performance anytime
- Ensure accurate assay execution
- Automate pipette checks to maintain standards
- Digitally manage tasks for improved lab efficiency

Ingenix Suite GxP Compliance Module

- Track system and instrument events for 21 CFR Part 11 and EMA Annex 11 compliance
- Access a comprehensive audit trail anytime
- Centralize audit trail data in an SQL database
- Digitally review and approve reports
- Review and approve reports digitally

Ingenix Suite Balance External Calibration Module

- Regularly check balance calibrations and manage weight sets digitally
- Verify calibration status per SOP
- Create bundles to track weight sets with balances
- Access all weight set information in one place

Ingenix Suite Advanced Pipette Calibration Module

- Digitally manage pipette calibrations for efficiency
- Calibrate per 2022 ISO8655 standards
- Use with 3rd party balances
- Automate tasks to boost efficiency

Further Highlights

- Automatic data transfer and storage in a secure SQL database
- Software detects data changes for integrity
- Authorization linked to user's Active Directory account
- Customizable role-based access control
- Available in 1-year and 5-year licenses

My Sartorius - Your Entry Point for Sartorius Digital Ecosystem

Your personalized dashboard will further provide you a tailored overview on all contents you may need. Interactive widgets allow you to access those Sartorius services relevant to you.

- Track your latest orders
- Access product certificates
- Download your purchased software
- Seamlessly transition to our eShop
- Media Center Exclusive for Partners
- Configure your new balance and request a quote
- Further apps to come soon

Comprehensive Solutions for Your Application Areas

Leverage Sartorius expertise built on a foundation of long-standing knowledge and continual innovation. With customized solutions for your application area, we provide everything you need for success in the Life Science, Biopharmaceutical Manufacturing, Quality Control and Testing, and Applied Industries.

Move your work forward with our portfolio and exceed your goals for reliability, reproducibility, and rapid results!

Our 2025 Spotlight Campaigns



Quality Control

Sartorius specializes in delivering innovative QC instruments and consumables for microbiological, physicochemical, and biological analyses, supporting accurate and reliable results. With a focus on efficiency and compliance, Sartorius empowers laboratories worldwide to streamline their Quality Control (QC) processes and maintain the highest standards in product safety and potency.



Biologics Discovery and Development

Biologic drugs have revolutionized medicine, offering hope for patients with serious conditions. Sartorius aids scientists by providing innovative solutions to enhance preclinical trials, boosting productivity and reducing drug development time and costs. Our products streamline purification processes, including cell culture clarification, antibody capture, polishing, concentration, and sterile filtration.

Sales and Service Contacts

For further contacts, visit
sartorius.com

Germany

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

USA

Sartorius Center of Excellence USA
3874 Research Park Drive
Tech Loop at the Ann Arbor Research Park
Ann Arbor, MI 48108, USA
Phone +173 4769 1600

If you have any questions, please contact your Sales
Representative or any of our authorized Sales Partners.

