



Highlights of
Our Lab
Solutions

Simplifying Progress

SARTORIUS

The Latest Solution 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.



Lab Weighing



Bioanalytics



Lab Filtration



Lab Water



Microbiological Analysis



Liquid Handling



Service & Training



Bioanalytics

Intellicyt® iQue3

Faster, Smarter Flow Cytometry

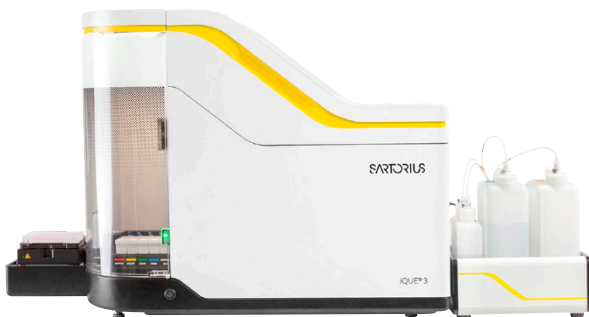
Perform rapid, multiplexed analysis of cells and beads in suspension with iQue3. Generate high-content data from small samples – ideal for screens where cells are precious and often very limited in number.

Features

- All-in-one ForeCyt® software workflow for data analysis and visualization
- Application-specific cell and bead-based assay kits
- Multiple instrument configurations to meet technical and budget requirements

Benefits

- Fast – processes 96 well plates in < 5 minutes and 384 well plates in < 20 minutes
- Data-rich results – multiplexed analysis of cells and beads in suspension for deeper insight
- Miniaturized assays – conserves precious cells and saves reagent costs



IncuCyte® Live-Cell Imaging

Explore Real-Time Automated Live-Cell Imaging and Analysis Inside Your Incubator

Derive deeper and more physiologically relevant information about your cells.

Features

- Real-time live-cell imaging and analysis
- Accommodates up to six microplates at a time – multiple users
- CMOS camera and efficient optics

Benefits

- View and analyze cells in real time
- Capture images over days and weeks
- Track experimental conditions from setup through analysis





Bioanalytics

Octet HTX

Select Antibody Lead Candidates with Confidence

Use the real time label-free Octet instruments to rapidly screen and select your antibody lead candidates with confidence while cutting down on assay time. Rapidly identify clones with high affinity and slow off-rates and monitor antibody titer in minutes and not hours.

Features

- Micro-fluidics free Dip and Read – reduces assay time and maintenance costs
- Flexible read heads – simultaneously reads 8, 16, 32, 48 & 96-wells
- Full plate off-rate rankings in minutes, not hours
- Automation ready

Benefits

- Compatible with crude supernatants – saves time and costs
- High-throughput – Analyze up to 96 antibody clones simultaneously for titer, off-rates and other CQAs
- Versatile – One instrument; multiple applications



Lab Water

Arium® Mini & Pro

Tailored to Your Specific Requirements

Arium® water system best suited for your sensitive analytical application.

Features

- Reliable ASTM Type 1 water quality
- Consistently low TOC levels
- Reliable process stability

Benefits

- Increase sensitivity of analytical results and minimize chemical ion interference
- Remove TOC levels down to a minimum and avoid background noises in chromatograms
- Offering a variety of services, e.g. installation, equipment qualification (IQ/OQ), preventative maintenance etc.





Liquid Handling

Picus® and Tacta® Pipettes

Tools for Reliable and Easy Pipetting

Sartorius simplifies pipetting and offers repeatable results time after time. The pipette and tips are designed to fit together and let you produce repeatable results at the level you need them to be.

Features

- Intuitive and simple user experience
- Low pipetting and tip ejection forces that reduce the risk of repetitive strain injury (RSI)
- Uniquely lightweight and compact
- Advanced safety features: Optilock, Safe-cone filters, password protection, calibration reminders

Benefits

- Achieve reliable results
- Be kind to your hand, by choosing Tacta and Picus pipettes
- Speedup your work with electronic Picus & Picus Nxt



Pipette Tips

Tips Designed to Match Your Pipette

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





Lab Weighing

Entris® II

Best Value in Its Class for Basic Weighing Tasks

No matter what you're weighing, the new Entris® II balance is always the right choice. Offering unrivalled value and backed by almost 150 years of German engineering expertise, the Entris® II comes in two product lines, so you can find the balance that meets your specific weighing needs.

Entris® II Essential line

Features

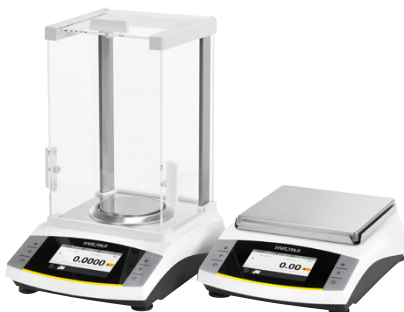
- isoCAL
- LED touch technology
- 12 built-in applications at a budget price.



Entris® II Advanced line

Features

- Real time level support
- Integrated protection systems
- CalAuditTrail
- Graphic touch display
- 13 built-in application programs



Cubis® II Balances

Data Integrity and Connectivity

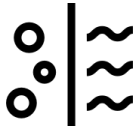
Need to comply with 21CFR11? Need the fastest stability on the market? Cubis® II – Your Balance, YourWay.

Features

- 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

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





Lab Filtration

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



Sartorius Minisart
Selection Guide:



Claristep® Filtration System

Sample Preparation of Eight Samples in One Go!

Do you process dozens of samples each day? Need to reduce the time, effort and waste in sample preparation? Claristep® - Ideal solution for sample preparation by particle removal prior to analytics.

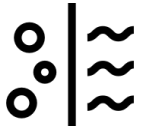
Features

- Multiplexing of up to 8 samples
- Syringeless filters and simple manual processing
- Filter sample volumes from 60 to 600 µl
- Made of highest purity materials

Benefits

- Fast and easy to use to maximize your throughput
- Reduced stress and waste with syringeless filtration
- Short sample contact time and high purity to achieve best analytical results





Lab Filtration

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 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.



Vivaspin® and Vivaflow®

Fastest Sample Concentration With the Highest Recoveries

The Vivaspin® and Vivaflow® ultrafiltration devices provide the most significant range of high recovery and super-fast protein concentrators for research and development labs. With the widest range of membrane and device options for all applications, including virus, mAb and exosome preparation.

Features

- 100µL to 5L concentration
- Diafiltration operation
- RC, PES and Hydrosart® membrane options

Benefits

- Highest recoveries with optimal membrane choice
- Fastest mL/min concentration available
- Reliable technical and application support





Microbiological Analysis

Rapid Microbiology Testing Kits

Better Be on the Safe Side Fast!

Bacteria, Fungi & Mycoplasma kits deliver results in only 3 hours. It is the perfect solution for the release of ATMPs, biopharmaceutical in-process testing or research purposes.

Features and Benefits

- PCR technique delivers results in only 3 hours
- Internal control reduces the risk of false positives
- TaqMan (R) probes provide high reaction specificity



Sterisart Sterility Testing 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





Microbiological Analysis

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 Air Monitoring Systems

Fixed or Portable Viable Air Monitoring for Controlled Environments

The MD8 Viable Air Monitoring systems combined with the gelatin membrane filters constitutes a sampling system for accurate, reproducible and quantitative detection of microorganisms & viruses. The samplers can also be used with BACTair culture plates for impaction sampling.

Features and Benefits

- Gelatin membrane filters are capable of detecting the smallest airborne viable contaminants, including viruses.
- Sampling head compatibility with in-line sterilization processes allows for disinfection and prevention of contamination.
- Continuous 8 hours sampling facilitates EU GMP Annex 1 regulatory compliance.





Service & Training

We Ensure the Quality of Your Results

At Sartorius, high-quality products go hand in hand with professional service. With our service offerings, we ensure the safe and optimal operation of your laboratory instruments. That means we cover the entire life cycle of your equipment from commissioning to qualification to regular maintenance.



Training Courses and Seminars

Sartorius offers a wide range of training courses that teach the theoretical fundamentals about Sartorius products and specific applications. They are designed to streamline your workflows, making you more efficient, while raising quality awareness.

For more information visit www.sartorius.com/service

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