

# Highlights of Our Lab Solutions

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.



Lab Weighing



Bioanalytics



Lab Filtration



Lab Water



Microbiological Analysis



Liquid Handling



Service & Training



## Bioanalytics

### iQue® High-Throughput Screening (HTS) by Cytometry

Truly Fast.  
Simply Effortless.

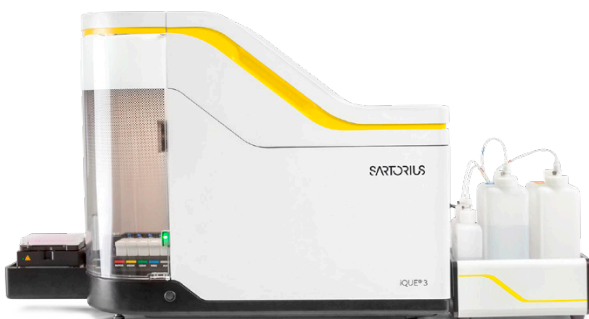
Perform rapid, multiplexed analysis of cells and beads in suspension with iQue®. Generate high-content data from small samples – ideal for screens where cells are precious and often very limited in number.

#### 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 Analysis

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® Label-Free Platforms

#### Best Performance, Exceptional Value

Don't sacrifice quality and performance when buying a label-free instrument for biomolecular interaction analysis. Sartorius is the only company to offer the choice of both industry-leading technologies – BLI and SPR. The entire Octet® range has been designed with one clear focus – to allow researchers to generate the most accurate data possible in the shortest timeframe imaginable.

##### Octet® BLI Features

- **High-throughput** – Analyze up to 96 samples simultaneously for binding kinetics, titer, off-rates and other CQAs
- **Crude Sample Compatibility** – Minimize time and costs analyzing unpurified samples
- **Low Total Cost of Ownership** – Low-maintenance fluidic-free system

##### Octet® SPR Features

- **Superior Gradient Injections** – Accurate measurement of molecule kinetics and affinity from a single analyte concentration
- **Optimized Fluidic System** – Minimized potential sources of blockage and benefit from efficient decontamination protocols
- **Independently Functioning Buffer Lines** – Use two running buffers during an assay



## Cell Selection and Retrieval

### CellCelector Flex

#### Automated Cell Selection and Retrieval System

With CellCelector Flex, you have everything you need for imaging, screening, picking and verifying target cells in one place. It's a versatile and automated all-in-one solution that reduces steps, streamlines workflows, and accelerates processing – while capturing cells without damage and enabling incredibly high viability.

##### Features and Benefits

- Picking and isolation from samples of single cells to complex models like organoids from a variety of liquid and semi-solid media
- Isolation of living or fixed cells
- Validated protocols for a wide range of applications
- Flexibility to develop unique imaging, detection and picking protocols
- Fully automated picking modes with real-time visualization of picking process
- Automated adjustment of picking position and height
- Complete documentation for traceability





## Liquid Handling

### Picus® 2 and Tacta® Pipettes

#### Tools for Reliable and Easy Pipetting

Sartorius simplifies pipetting and offers repeatable results time after time. The Sartorius pipettes and tips are designed to a complete system and let you produce repeatable results at the level you need them to be. Experience perfect pipetting utilizing superior ergonomics, efficiency and safety.

##### 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® 2 pipettes
- Speedup your work with electronic Picus® 2



### Pipette Tips

#### Tips Designed to Match Your Pipette

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

##### Features

- Perfect and soft sealing with Sartorius pipettes
- Unique safety space between filter and sample
- Packing options for every lab and color-coding to match your pipette

##### Benefits

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





## Lab Weighing

### Quintix® Pro

#### Redefining the Standard

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

- Ensure the highest weighing accuracy with the internal motorized levelling (optional) and automatic internal adjustment isoCAL
- Easy access to menu option with a large, 7-inch, high resolution full-touch graphic display with modern upgraded user interface
- Smart design and chemically-resistant surfaces made of PBT, glass and stainless steel support routine cleaning protocols
- User-defined, automatic start timer powers up the instrument in time for daily use
- Quick access to the integrated digital manual for common help topic
- 15 integrated application programs for common weighing tasks



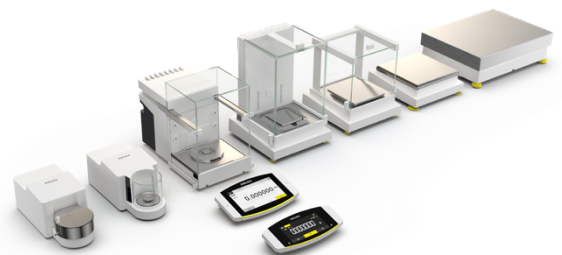
### Cubis® II Balances

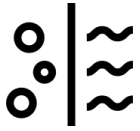
#### Data Integrity and Connectivity

Weighing in a regulated environment can be complex, with stringent compliance directives and controls required to deliver the accuracy and reliability of results required by the pharmaceutical industry. One of the technologies helping scientists to achieve this level of accuracy is the Cubis® II balance series from Sartorius.

##### Features with display unit MCA

- Thousands of hardware and software configurations supporting your personal preferences
- Seamless connectivity and digital data management to diverse laboratory information systems without the need of additional software
- Full end-to-end data integrity. 21 CFR Part 11 compliance, integrated audit-trail, state-of-the-art user management and electronic signature
- Full data integrity following ALCOA+ principles
- User management including centralized password management support ("single-sign-on" via LDAP), high security password configuration, login rules
- Software application with guided workflow for routine and deep cleaning
- Outstanding service support. Integrated status center, integrated service functions preventive maintenance on accredited standards





## 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 cm<sup>2</sup>
- 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:



### Membrane Filters

#### High Efficiency Membranes for Filtration and Venting Applications

Sartorius produces a wide variety of microporous membranes that are especially designed, developed and manufactured to meet differing needs of the biotechnology industry.

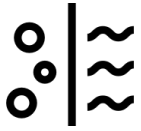
##### Features

- High flow rates
- Broad and strong chemical compatibility
- Low content of extractables

##### Benefits

- Sartorius core competency
- Covers essential applications in the life science industry
- Highly individual product range
- Available in a wide range of diameters and membrane types





## 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 1 L 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.



### Vivaspin® and Vivaflow®

#### Efficient Concentration and Buffer Exchange of Any Molecule

Vivaspin® and Vivaflow® represent the most comprehensive range of ultrafilters for sample concentration and buffer exchange. Offering state-of-the-art designs and a wide selection of ultrafiltration membranes, these products provide unparalleled speed and flexibility to reliably process any protein, nucleic acid, virus, nanoparticle or other macromolecule.

##### Features

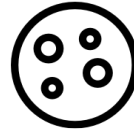
- TFF technology, advanced flow channel designs, and intuitive handling
- Multiple membrane chemistries and feed capacities from 0.1 mL to 5 L
- Low-volume, pipette-friendly dead stop pockets
- Single use solutions that need no additional cleaning steps

##### Benefits

- Reduce processing time, eliminate blocking and forget process optimization
- Ensure maximum recovery of any molecule from any sample volume
- Prevent sample losses while achieving concentration factors up to 100X or more
- Maintain sample integrity and improve laboratory safety







## Microbiological Analysis

## Sartobind® Lab & Vivapure®

### Faster and More Flexible Biomolecule Purification

Sartobind® membranes are available in a choice of consumables and kits that ensure flexible and reliable biomolecule purification with or without equipment. By switching to membrane chromatography, you can eliminate many of the limitations of resin columns, to condense workflows and significantly reduce your time-to-result.

#### Features

- Plug and play units are ready-for-use
- Spin and pressure filter formats suit multiple handling methods
- Macroporous membranes with high-density IEX and affinity ligands
- Process scale capsules and cassettes available with the same membranes

#### Benefits

- Eliminate column packing and buffer degassing
- Reliable purification with or without equipment
- Purify higher quantities of antibodies, other proteins or viruses in seconds, not minutes
- Seamless transition from research to clinical and commercial production



## Rapid Microbiology Testing

### 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® probes provide high reaction specificity





## Microbiological Analysis

### Sterisart® Sterility Testing System

#### Where Convenience Meets Compliance

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
- Patented dual needle design
- Go fully paperless with the Sterisart® Gen 4
- Built for 21 CFR Part 11
- Semi-automation & visual guidance of workflows



### 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.





## Lab Water

### 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.



### 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.





## Service & Training

# We Ensure the Quality of Your Results

### Your Success Is Our Mission

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 on Sartorius Services please visit:  
[www.sartorius.com/services](http://www.sartorius.com/services)



#### 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

 **Find out more:** [sartorius.com](http://sartorius.com)