

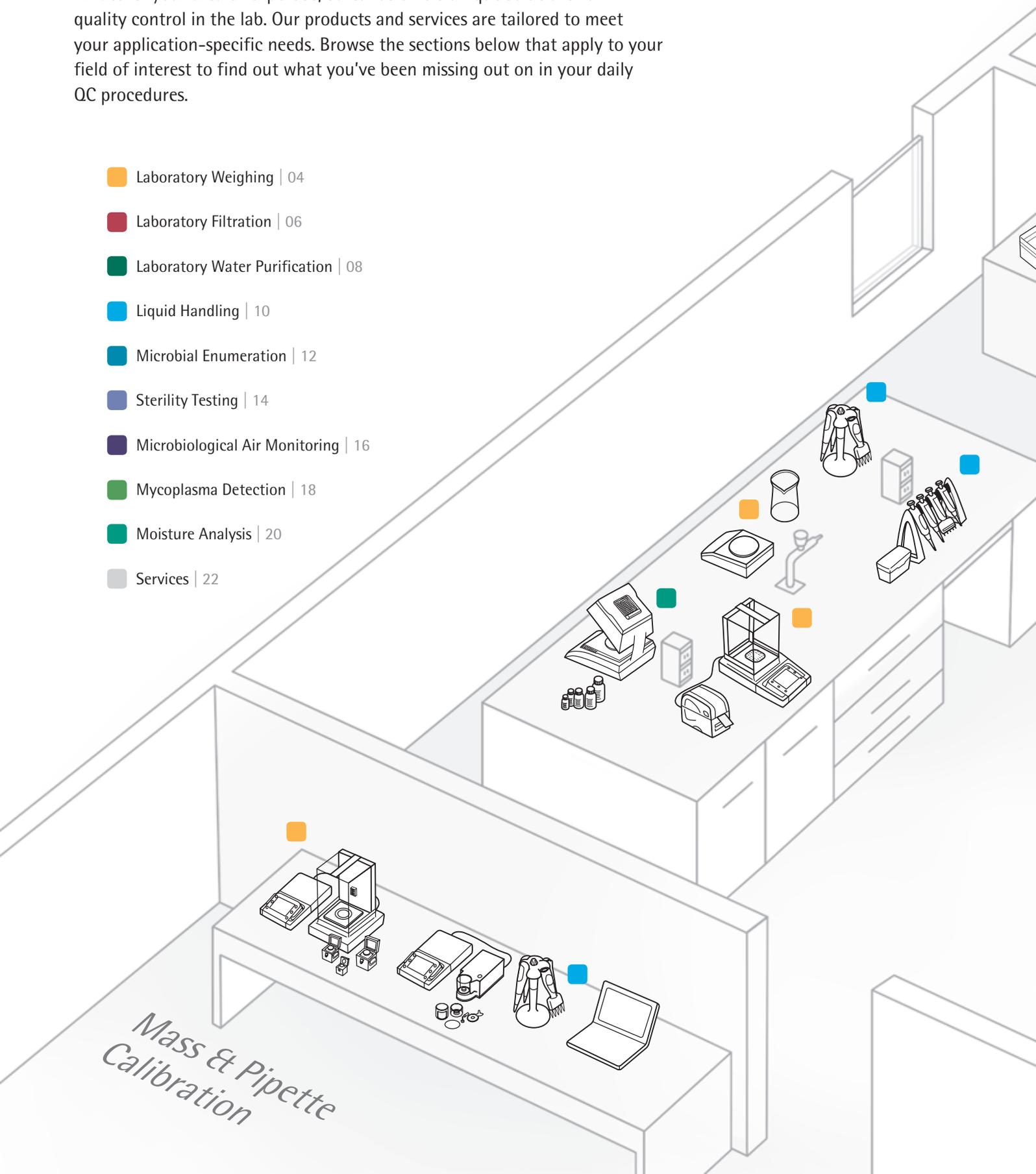
## Quality Control in Pharmaceutical Laboratories Our Product Solutions for Your Needs



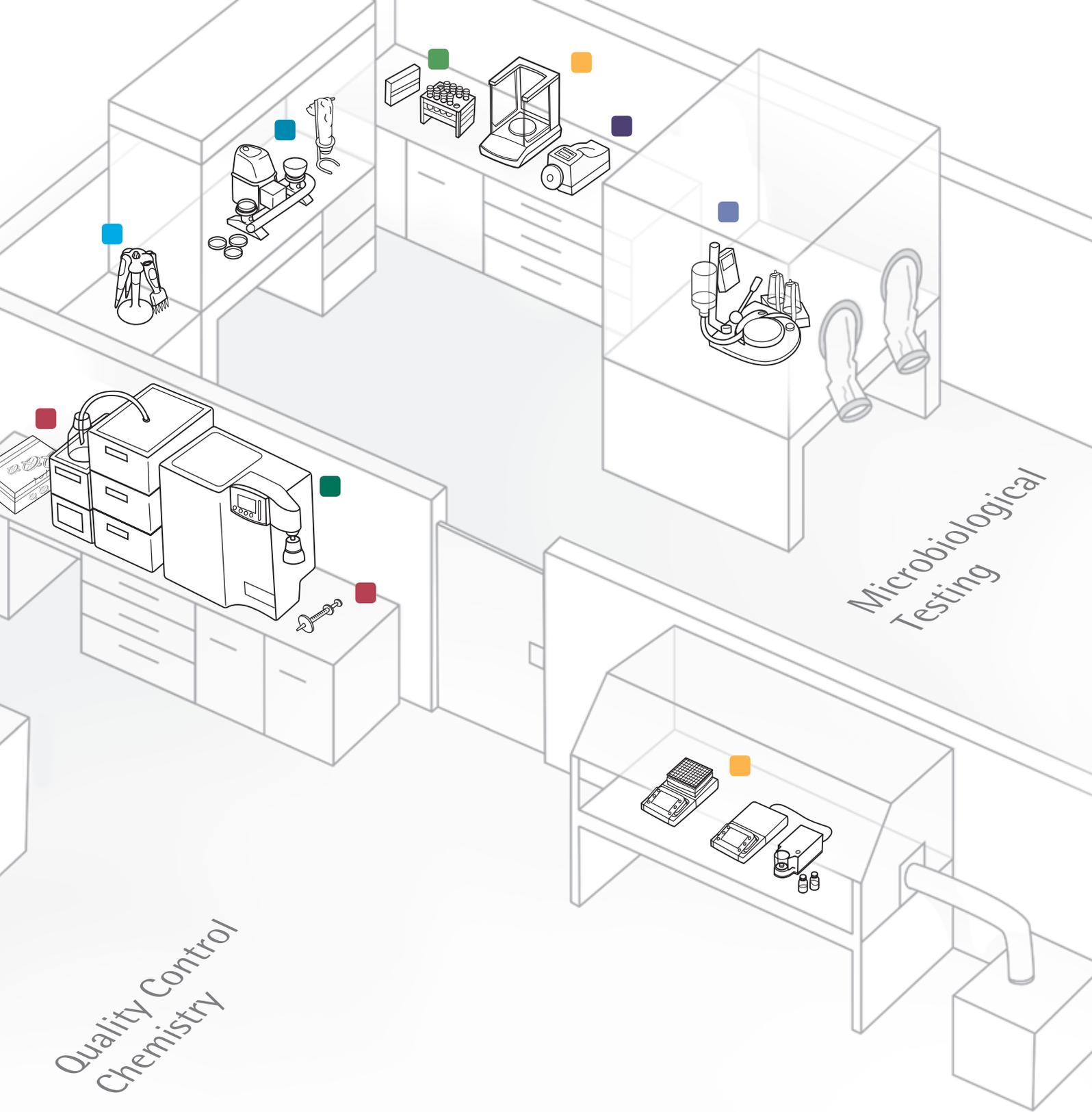
# Discover the Potential of Your QC Lab

Whatever your area of expertise, Sartorius offers unique solutions for quality control in the lab. Our products and services are tailored to meet your application-specific needs. Browse the sections below that apply to your field of interest to find out what you've been missing out on in your daily QC procedures.

-  Laboratory Weighing | 04
-  Laboratory Filtration | 06
-  Laboratory Water Purification | 08
-  Liquid Handling | 10
-  Microbial Enumeration | 12
-  Sterility Testing | 14
-  Microbiological Air Monitoring | 16
-  Mycoplasma Detection | 18
-  Moisture Analysis | 20
-  Services | 22



Mass & Pipette  
Calibration



Quality Control  
Chemistry

Microbiological  
Testing

### Services for Your Laboratory

- Installation
- Extended warranty
- GMP | GLP qualification (IQ | OQ)
- Calibration services
- Preventive maintenance
- Service contracts
- User training | Laboratory Academies

## ■ Laboratory Weighing

Configure your balance to suit your individual needs and seamlessly integrate it into your process. Our modular-designed series of Cubis® lab balances enables you to combine your favorite display and control unit with a weighing module, draft shield, data interface and many more features of your choice.

### Speed

- Obtain lightning-fast display response to accelerate your weighing procedures.
- Speed up your system setup and calibration.
- Benefit from flexible data acquisition and communication capabilities.
- Reduce your validation time thanks to complete documentation and support.

### Regulatory Compliance

- Meet the requirements of ISO, GMP, GLP, FDA 21 CFR Part 11, USP and other regulations.
- Obtain completely traceable results according to the recommendations of the International Organization of Legal Metrology (e.g., OIML R76-1) and national standards.
- Use the Advanced Pharma Compliance (APC) balance functionality for reliable process monitoring and qualified results.

### Ease of Use

- Easily operate your balance with interactive user guidance and adapt it to your needs.
- Configure tasks for different users with a user | task management tool for consistent, error-free results.
- Readily adapt the balance for your workflows.

### Quality of Results

- Use a balance that exactly meets your individual process, accuracy and quality requirements.
- Benefit from the highest level of weighing performance and technology.
- Reduce the influence of ambient conditions and different users with the help of self-tests and warning functions.



## Process Integration

- Easily integrate your balance into your lab processes and IT infrastructure with its flexible communication concept.
- Customize your balance with downloadable Q-Apps.
- Implement your SOPs using customer-specific Q-Apps.
- Take advantage of advanced Web Services communication with your LIMS, ELN, etc.
- Work with smart accessories for special applications.

## Applications

- Sample preparation, e.g. for HPLC, elemental analysis, thermal analysis
- Backweighing and filter weighing
- General weighing and density determination
- Compounding, totalizing, formulation
- Tablet checking
- Animal weighing
- In-process control | PC
- Calibration of weights
- Pipette checking and calibration

## Products



**Cubis® with an MSA Display**  
Offers the highest level of convenience. Provides optimal support in regulated areas. Integrates into your customized process workflows with downloadable Q-Apps.



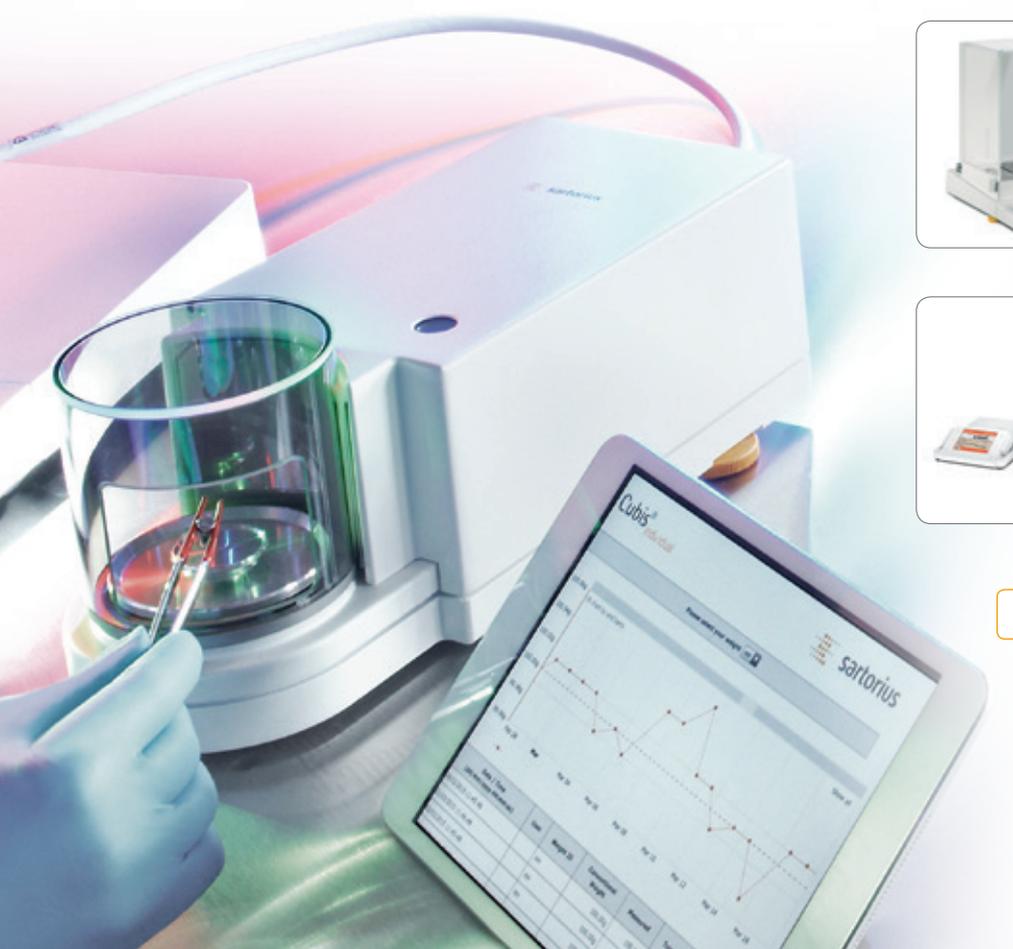
**Cubis® with an MSE Display**  
Perfect for focusing on the essentials of weighing: high-precision weighing with logical and straightforward operation.



**Cubis® MCM Manual Mass Comparators**  
Cover capacities from 6.1 g to 64 kg and readabilities from 0.1 µg to 10 mg. All models have climate sensors for temperature, humidity and air pressure and integrated workflow control.

## Services

- Installation
- Extended warranty
- GMP|GLP qualification (IQ|OQ)
- Calibration service (USP, ISO 17025)
- Preventive maintenance
- Service contracts
- User training | seminars



## ■ Laboratory Filtration

Get the exact type of syringe filter you need. Our Minisart® filters are ideal for a wide range of applications. They are a clean, safe and reliable choice for all of your low-volume and medium-volume filtration steps.

### Ease of Use

- Choose the right filter and minimize the loss of your samples. Ready-to-use Minisart® syringe filters for analytics are available in three different sizes for sample volumes of less than 100 µL to 100 mL.
- Use filtration for the fastest liquid sterilization. It removes all microorganisms and particles reliably, maintaining the composition of the sample thanks to exceptionally low non-specific binding.

### Quality of Results

- Choose optimized filters for your sample preparation prior to analytics. Their polypropylene (PP) housing withstands harsh solvents and chemicals without adding leachables to the samples.
- Utilize filters made of medical-grade acrylic (MBS) for the sterile filtration of drugs and other aqueous solutions.
- Visit our Minisart® production site: All components are manufactured and quality controlled by Sartorius. Audits can also be ordered and performed by a Notified Body.



## Regulatory Compliance

- Meet the current GMP standards with Minisart®: Our quality management system has been certified for compliance with DIN EN ISO 9001 and DIN EN ISO 13485.
- Perform biosafe filtration according to USP Class VI Plastic Tests. In addition, comply with Title 21 of the Code of Federal Regulations, Sections 210.3(b)(6) and 211.72, for non-fiber-releasing filters. Minisart® NML is also classified in class IIa of the Medical Devices Directive 93 | 42 | EEC.

## Applications

- Sample preparation prior to analytics
- Clarification of liquids prior to downstream tests
- Sterilization of liquids prior to downstream tests

## Products



### Minisart® RC 25

Contains a regenerated cellulose (RC) membrane ensuring the lowest non-specific binding.



### Minisart® SRP 15

Has a PTFE membrane for the highest chemical compatibility and a pH range from 1 to 14.



### Minisart® NML

The surfactant-free cellulose acetate membrane is perfect for aqueous liquids.

## Services

- Validation guide
- CONFIDENCE® Validation Services including extractables | leachables studies, physiochemical studies and microbiological studies
- Lab Filtration Academy



# ■ Laboratory Water Purification

Find the perfect lab water system to meet all of your requirements.

To ensure you obtain the water quality you need, the new arium<sup>®</sup> product family gives you a choice of 70 system variations, each tailored to a specific application. Thanks to their modular design, arium<sup>®</sup> systems can be integrated into any lab environment.

## Regulatory Compliance

- Meet and exceed the ASTM Type 1 water quality standards.
- Fulfill pharmacopeia requirements for pure and ultrapure water.
- Rely on lab water systems manufactured in compliance with ISO 9001.

## Ease of Use

- Easily navigate using the touch-activated, high-quality glass display – even while wearing gloves.
- Monitor all important parameters at once utilizing the clearly organized user interface and the intuitive menu.
- Optimize your purified water quality with the intelligent iJust software. It improves your recovery rate for better system performance and economical use of resources.
- Keep your water usage safe and reliable thanks to visual alert management.
- Eliminate time-consuming and labor-intensive cleaning procedures by using disposable bags instead.

## Quality of Results

- Obtain reproducible and reliable results in analytical applications by efficiently reducing TOC to less than 2 ppb.
- Remove endotoxins, microorganisms, DNases and RNases safely and permanently, based on ultrafiltration by the recirculation loop.
- Count on constant water quality of the highest level due to top-down flow cartridge technology and carefully selected ion-exchange resin.
- Rely on safe system operation supported by highly accurate TOC monitoring.
- Safely store your water in single-use bagtanks made of pharmaceutical-grade materials.



## Flexibility

- Save space thanks to the compact design: Adapt your water system to any lab environment by flexibly positioning the display and dispensing unit.
- Maximize available work space with an easy-to-use remote dispenser that is designed to fit all benchtop, wall-mounted and built-in units.
- Conveniently move your stored purified water with the flexible bagtank on rollers.
- Custom-tailor your system on a modular basis for a perfect match designed for your specific application.

## Applications

- Analytical Methods:  
HPLC, GC-MS, AAS, ICP-MS, IC, immunocytochemistry, endotoxin analysis, TOC analysis, photometry, ELISA, RIA
- Preparative Methods:  
Buffer preparation, PCR, electrophoresis, photometry, media for cell cultures, production of monoclonal antibodies, histology sections
- Feed Lab Instruments:  
Autoclaves, labware dishwashers, humidifiers, water baths, rinsing vessels, etc.

## Products



**arium® Lab Water Systems**  
The arium® line offers a wide range of modular-designed systems for producing ASTM Type 1 ultrapure and pure water.



**arium® bagtank**  
Stores pure water in a single-use bag. The bagtank reliably protects purified water from secondary contamination. It is the ideal solution for time-saving and reliable operation.

## Services

- Installation
- Extended warranty
- GMP|GLP qualification (IQ|OQ)
- Exchange of consumables
- Calibration service
- Preventive maintenance
- Service contracts
- User training | seminars



## ■ Liquid Handling

Obtain reliable pipetting results effortlessly and enjoy your work. Our Picus® and Picus® NxT electronic pipettes, as well as the mechanical mLINE®, offer excellent usability, exceptional ergonomics and high reliability of results – especially when used with Sartorius pipette tips.

### Regulatory Compliance

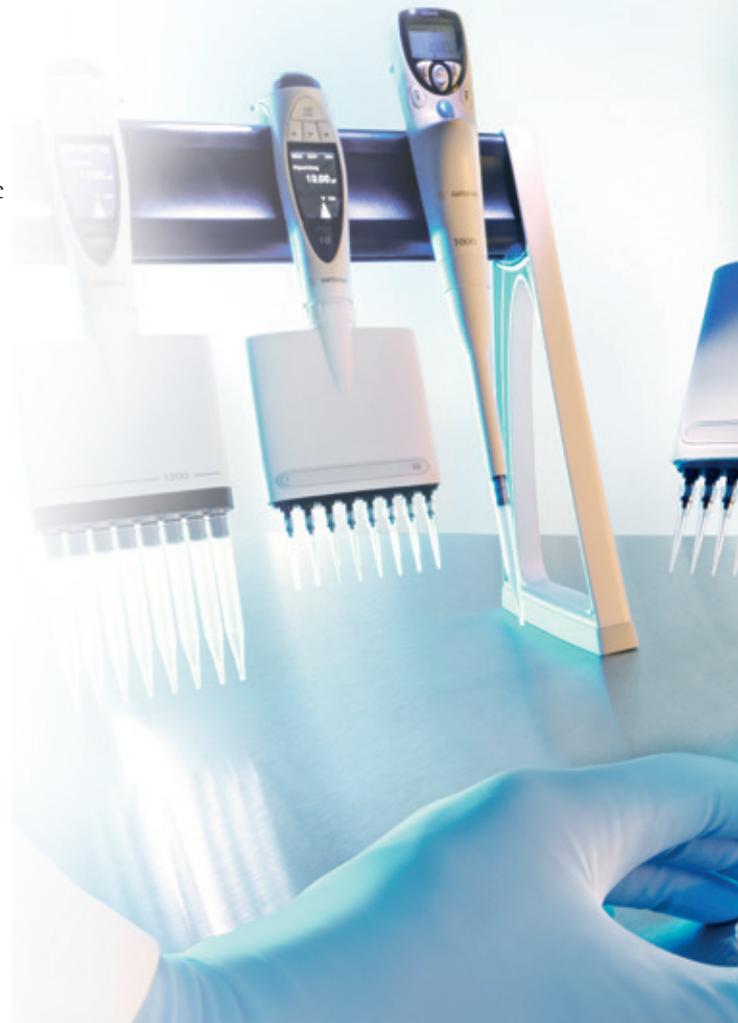
- Comply with ISO 8655 by using our liquid handling products and calibration services.
- Receive lot-specific purity certificates for tips.
- Conform to the strictest lab regulations with the Picus® NxT certificate of accredited 3-point calibration.
- Rely on our calibration services with ISO 17025 accreditation.

### Ease of Use

- View the volume quickly and easily on the large, crisp display.
- Intuitively navigate the clear user interface of our electronic pipettes.
- Make pipetting tasks quick and easy by using the extensive range of pipetting modes offered by Picus® and Picus® NxT.
- Quickly clean and maintain your pipettes designed for easy disassembly for servicing.
- Easily find the matching color-coded tips and pipettes.

### Quality of Results

- Use Sartorius pipettes and tips for the best accuracy and precision.
- Easily calibrate and adjust your pipettes for optimal results with various liquids.
- Prevent contamination with cost-effective tip-cone filters.
- Obtain reliable results for liquids with low surface tension by using low retention tips.



## Ergonomics

- Work effortlessly and ergonomically – no matter whether you are left- or right-handed.
- Prevent repetitive strain injury (RSI) with the help of the ergonomic design and low pipetting forces of our pipettes.
- Pick up the tips with no effort thanks to the Optiload feature with spring-loaded tip cones.
- Eject tips easily with the electronic tip ejection of Picus® and Picus® NxT.

## Applications

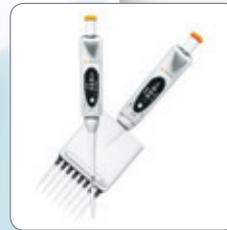
- Analytical Methods:  
HPLC, immunochemistry, endotoxin analysis, photometry, LAL test, ELISA, etc.
- Preparative Methods:  
Buffer preparation, PCR methods, electrophoresis and media for cell cultures, etc.
- General Liquid Handling:  
Routine pipetting, dispensing, sample and reagent transfer, dilution, etc.

## Products



### Picus® and Picus® NxT

The smallest and lightest electronic pipettes on the market are comfortable to hold and speed up your work with their wide range of pipetting modes.



### mLINE®

A mechanical pipette that provides excellent ergonomics and thus maintains high accuracy and precision also during long pipetting series.



### Pipette Tips

Designed to match Sartorius pipettes for the best possible results. Choose from standard, filter or low retention tips in racks, refill towers, packs or packages.

## Services

- ISO 8655 calibration
- ISO 17025 calibration certificates
- Preventive maintenance | repairs
- Pipetting Academy:  
Training in ergonomics, pipetting techniques, calibration and servicing of pipettes



## ■ Microbial Enumeration

Simplify your microbiological quality control. Our Microsart® consumables and hardware products enable you to streamline your procedures and to work in an ergonomic position. Their accurate and dependable test results help you meet the challenges and demands of today's healthcare environment.

### Regulatory Compliance

- Perform microbial limit tests using our Microsart® @media media dishes to ensure 100% compliance with EP 2.6.12 and USP <61>.
- Obtain certification on sterility, growth promotion, retention as well as chemical and physical parameters, including pH and flow rate.

### Ease of Use

- Remove the funnel fast and easily after filtration thanks to its practical click-fit closure.
- Monitor the progress of your filtration run just by viewing the transparent funnel.
- Effortlessly transfer your membrane filter onto the agar dish and mitigate the risk of RSI.
- Streamline your workflows and simplify testing with ergonomically friendly consumables and hardware.

### Quality of Results

- Eliminate membrane manipulation by touch-free membrane transfer. This will minimize a major source of secondary contamination.
- Prevent secondary contamination during assembly with a single-use filtration unit that is sterile and ready to use.
- Thoroughly rinse your funnel after filtration without retaining any liquid due to its optimized design.



## Speed

- Save time with only a few simple steps from filtration to incubation.
- Forget about the hassle of assembly with our ready-to-use filtration unit combining funnel, base support and filter.
- Transfer the membrane in a sterile manner without tweezers.
- Improve your time to results.

## Applications

Microbiological quality control for:

- Water testing
- Bioburden testing
- Raw material testing

## Products



### Microsart® @filter

Ready-to-use sterile filtration unit that combines a funnel, a filter base and a gridded membrane filter. Various pore sizes and filter colors for contrasting backgrounds are available.



### Microsart® @media

Prefilled media dishes that allow touch-free membrane transfer– without the need for tweezers. This reduces the risk of false-positive results.



### Microsart® e.jet

Vacuum pump with a diaphragm membrane. It is designed to create sufficient vacuum for filtration and to allow simultaneous transfer of the filtered liquid to waste.

## Services

- GMP|GLP qualification (IQ|OQ)
- Preventive maintenance



## ■ Sterility Testing

Obtain convincing proof of the performance of our innovative solutions for sterility testing of pharmaceuticals using membrane filtration. Our Sterisart products fully comply with international pharmacopeias.

### Regulatory Compliance

- Get Sterisart disposables for your sterility tests that are fully compliant with international pharmacopeias, such as USP <71>, Eu. Ph. 2.6.1 and JP 4.06.
- Receive full documentation, including IQ/OQ and validation guides.
- Ensure GMP compliance with quality assurance certificates and individually traceable units.
- Comply with ISO 14644 thanks to our particle-emission-free Sterisart pump.

### Ease of Use

- Conduct easy and reliable aseptic sampling with our unique Septum version – without any risk of contamination or accidental injuries. You no longer need to cut the tubes or pierce the tubing with a syringe.
- Easily find and grasp the large, 4-edged red caps – even while you are wearing a double pair of gloves. The caps are conveniently connected to the barb of the sterilizing-grade vent filter.
- Quickly clean and maintain your Sterisart NF Universal pump due to its easy-to-disassemble design.
- Rely on process traceability via barcode scanner: Installed on an upgraded pump version, it provides convenient, error-free data transfer and storage (batch and individual number).

### Quality of Results

- Minimize the risk of operating errors with color-coded tubing clamps and a barcode on the packaging.
- Sample, inject and dilute with medical-grade septum versions to avoid taking risks.
- Benefit from the highest accuracy with proven equal distribution of liquid to both containers.



## Speed

- Install Sterisart tubing fast and easily in your Sterisart pump – without having to interrupt the pumping process.
- Run the highest pump speeds without excessive foaming due to the inlet design of the test containers.
- Save time by using the dual-needle spike with a sterilizing-grade filter instead of a separate venting needle.

## Applications

- Routine sterility testing in cleanrooms and isolators
- Wide variety of sample types: liquids, powders and emulsions
- Sterility testing of a broad range of sample vessels: bags, syringes, bottles, vials and blow-fill-seal devices

## Products



### Sterisart NF

Choose from more than 20 different single-use filtration units. Test the sterility of liquids, powders and antibiotics in open or closed containers.



### Sterisart Universal Pump

This low-maintenance, H<sub>2</sub>O<sub>2</sub>-resistant AISI 316 L stainless steel pump equally distributes sample and media in the Sterisart containers for sterility testing. The high-end version guides you by individual menu prompts.

## Services

- Validation and revalidation (H<sub>2</sub>O<sub>2</sub> gas tightness; bacteriostasis | fungistasis testing)
- Installation
- Extended warranty
- GMP|GLP qualification (IQ|OQ)
- Preventive maintenance
- Service contracts
- User training | seminars



## ■ Microbiological Air Monitoring

Discover our field-proven air samplers. They are routinely used for active air sampling as well as the quantitative detection of airborne microorganisms and even viruses. With our matching ready-to-use membrane filters and media plates, you will obtain reliable results fast and easily.

### Speed

- Easily install the gelatin membrane filter and BactAir® impactation plates by simple attachment to the adapter.

### Ease of Use

- Forget about cleaning validation: All consumables are single use, including the sampling head.
- Benefit from ready-to-use disposables: single- and triple-packaged and gamma-sterilized consumables are available.
- Reduce the risk of secondary contamination by transfer of the gelatin membrane filter – without using tweezers – to the culture medium plate.
- Monitor controlled environments with our selection of portable, battery-powered as well as stationary air samplers or integrate MD8 airscan samplers into isolators, blow-fill-seal machines and filling lines.

### Quality of Results

- Achieve the highest recovery rates while sampling airborne microbes and particles with the gelatin filtration method.
- Safely sample yeasts, fungi, bacteria and even viruses.
- Sample small microbes and particles with the impactation method on agar plates thanks to the optimized design of the single-use sampling head.



## Regulatory Compliance

- Rely on full compliance with international regulations, e.g., USP 1116, EN ISO 14644 and EN ISO 14698, with the MD8 air sampler and materials for microbiological monitoring of airborne particles.
- Ensure GMP compliance with quality assurance certificates.
- Perform isokinetic air sampling required by (EU) GMP guidelines | EN ISO 14644 – without interfering with zone protection.
- Calibrate the air flow for accurate sampling of 1 m<sup>3</sup> of air as required by regulations.

## Applications

- Environmental air monitoring of unspecified zones, controlled zones and cleanrooms
- Environmental air monitoring of highly sensitive and critical applications in Grade A zones, such as isolators, filling lines and blow-fill-seal machines
- Washing out of antimicrobial agents, such as antibiotics and cytostatics

## Products



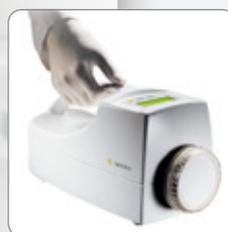
### Gelatin Membrane Filters (GMF)

The sampling head along with the gelatin membrane filter contained is presterilized and ready to use for differentiated detection of airborne microbes.



### BactAir® Culture Media Plates

Ready to connect as collecting heads to AirPort MD8, these prefilled and presterilized plates are available in a choice of media (total count or detection of fungi).



### AirPort MD8

Portable, battery-powered air sampler for use with gelatin membrane filters (GMF) and BactAir® culture media plates.

## Services

- GMP|GLP qualification (IQ|OQ)
- Preventive maintenance
- User training | seminars



## ■ Mycoplasma Detection

Detect mycoplasma contamination at the earliest stage possible.

Our Microsart® Mycoplasma Detection Kits enable you to reduce your test time to only a few hours and achieve absolutely reliable results. Depending on your application and required safety levels, you can choose between two different kits.

### Regulatory Compliance

- Rely on Microsart® AMP and ATMP Mycoplasma that are validated for sensitivity, specificity and robustness according to EP 2.6.7 in combination with EP 2.6.21.
- Use our validation planning and reporting services for easy implementation.

### Ease of Use

- Save time and money with just a few simple steps from DNA extraction to interpretation of PCR results.
- Identify the various tubes easily and conveniently by their colored caps.
- Utilize the kits' tested, ready-to-use master mix with all the essential components including polymerase. You won't need an expert at your side.
- Benefit from the highest flexibility with different protocols for various sample materials.

### Quality of Results

- Minimize the risk of false-positive results by using TaqMan® probes to add specificity to your PCR test system.\*
- Utilize an internal and a positive control to decrease the risk of false-negative results.
- Minimize the occurrence of false-positive results. The kits contain dUTP instead of dTTP, so the option is available to degrade amplicons from previous analysis by use of uracil-DNA glycosylase (UNG).

\* TaqMan® is a registered trademark of Roche Molecular System Inc.



## Speed

- Reduce your testing time from 28 days to 3 hours.
- Achieve optimal results thanks to the flexible volume range of Microsart® AMP Mycoplasma from 200 µL to 18 mL. If you need highest sensitivity, use our centrifugal concentrator Vivaspin® 6 or 20 spin columns for concentration.

## Applications

- For analyzing cell culture materials and raw materials, for in-process controls and for lot-release testing of finished products
- Mycoplasma detection in ATMPs (advanced therapy medicinal products).

## Products



**Microsart® AMP Mycoplasma**  
With its very flexible sample volume (up to 18 mL) and sample preparation, the kit is specifically designed for the quality control of raw materials, in-process control and lot-release testing.



**Microsart® ATMP Mycoplasma**  
Especially designed for fast and easy detection of mycoplasma in advanced therapy medicinal products, ATMPs.

## Services

- User training | seminars
- Validation support



## ■ Moisture Analysis

Choose a moisture analyzer that matches the specific requirements of your pharmaceutical process and quality control applications. From our basic model up to our high-end solutions, all of our moisture analyzers are especially user-friendly and deliver fast and accurate results.

### Regulatory Compliance

- Use ASTM-compliant methods.
- Meet the requirements of quality standards, such as ASTM.

### Ease of Use

- Easily operate your moisture analyzer by following the guidance prompts provided by the interactive menu.
- Benefit from the modular system for more flexibility and higher efficiency.
- Implement up to 300 individual methods.

### Quality of Results

- Achieve the highest level of accuracy and measurement speed.
- Obtain reliable results with individually optimized methods.
- Utilize a reliable method with guaranteed comparability to approved reference methods.



## Speed

- Increase your flexibility with fast measurements.
- Connect up to 4 heating units for higher throughput.
- Set it up quickly: Select your product list and start.

## Applications

- Moisture analysis of pharmaceutical ingredients
- Analysis of the solid content of solutions
- Determination of the moisture content of drug formulations
- General moisture determination, even of very dry samples

## Products



### Mark3 HP, Mark3 LTE

The highest precision and performance for demanding applications. Their modularity makes them absolutely efficient.



### MA160

Manages sophisticated tasks. Ideal for performing fast and accurate analyses of a diverse range of product samples.



### MA37

This fast, reliable and robust device is ideal for routine moisture determination.

## Services

- Method development
- Extended warranty
- GMP|GLP qualification (IQ|OQ)
- Calibration service
- Service contracts
- User training | seminars



## ■ Sartorius Services

Extend the lifetime of your equipment with our comprehensive range of services. In order to keep your pharmaceutical lab working successfully, accurately and efficiently, we provide all the means to ensure highest reliability and uptime of your equipment as well as confidence in your results.

### Trained Operators

- Have us instruct your operating staff as part of each installation and IQ/OQ.
- Gain theoretical knowledge and acquire the skills necessary to put it into practice: Our application specialists and training center offer academies and seminars that will show you how to work even more efficiently and confidently.

### Optimized System Performance

- Ensure full performance right from the start and for the entire equipment lifecycle with our professional system installation.
- Get the most out of your equipment with our individual configuration and on-site adaption.

### Regulatory Compliance

- Use our installation and operational qualification (IQ/OQ) services to ensure that your equipment can be used in a highly regulated environment.
- Obtain documented quality and traceability of your measurement results in compliance with ISO 17025, GLP|GMP and FDA regulations as part of our calibration services. Our ISO 17025-accredited certificates are accepted worldwide and enhance your recognition as a qualified operator.

### Quality of Results

- Trust the accuracy and preciseness of your analytical results with regular equipment calibration. Sartorius will perform and document this service for you. We offer a variety of accredited and ISO calibration certificates which meet the requirements of the pharmaceutical and regulated industry.



## Range of Services

### Trouble-free Operation

- Rest assured that your equipment will run efficiently throughout its maximized lifetime.
- Ensure optimal equipment performance and prevent unnecessary interruptions or malfunctioning with our preventive maintenance.
- Benefit from the highest repair quality and minimize equipment downtime thanks to our professional repair services with original spare parts.



#### Installation Services

Our factory-trained service personnel ensures full system performance right from the start.



#### Extended Warranty

Extended warranty packages protect your investment and provide full cost transparency.



#### System Qualification (IQ/OQ)

Comprehensive and product-specific documentation will provide proof that your equipment operates in compliance with ISO 17025, GLP|GMP and FDA regulations.



#### Preventive Maintenance and Service Contracts

Special service contracts for your confidence and instrument reliability will help you achieve the highest output, consistency and efficient instrument usage.



#### Calibration Services

Documented quality and traceability of measurements ensure that your results will be compliant with ISO 17025, GLP|GMP and FDA regulations.



#### User Training | Academies

A wide selection of extensive training programs and academies will boost the successful, reliable and efficient work of your staff.



# Sales and Service Contacts

For further contacts, visit [www.sartorius.com](http://www.sartorius.com)

## Europe

### Germany

Sartorius Lab Instruments  
GmbH & Co. KG  
Weender Landstrasse 94-108  
37075 Goettingen

Phone +49.551.308.0  
Fax +49.551.308.3289

### France & Suisse Romande

Sartorius France  
2, rue Antoine Laurent de Lavoisier  
ZA de la Gaudrée  
91410 Dourdan

Phone +33.1.70.62.50.00  
Fax +33.1.64.59.76.39

### Austria

Sartorius Austria GmbH  
Modecenterstrasse 22  
1030 Vienna

Phone +43.1.7965760.0  
Fax +43.1.7965760.24

### Belgium

Sartorius Belgium N.V.  
Rue Colonel Bourg 105  
1030 Bruxelles

Phone +32.2.756.06.90  
Fax +32.2.481.84.11

### Finland & Baltics

Sartorius Biohit Liquid Handling Oy  
Laippatie 1  
00880 Helsinki

Phone +358.9.755.951  
Fax +358.9.755.95.200

### Hungary

Sartorius Hungária Kft.  
Kagyló u. 5.  
2092 Budakeszi

Phone +3623.457.227  
Fax +3623.457.147

### Ireland

Sartorius Ireland Ltd.  
Unit 41, The Business Centre  
Stadium Business Park  
Ballycoolin Road  
Dublin 11

Phone +353.1.8089050  
Fax +353.1.8089388

### Italy

Sartorius Italy S.r.l.  
Viale A. Casati, 4  
20835 Muggiò (MB)

Phone +39.039.4659.1  
Fax +39.039.4659.88

### Netherlands

Sartorius Netherlands B.V.

Phone +31.30.60.53.001  
Fax +31.30.60.52.917

[info.netherlands@sartorius.com](mailto:info.netherlands@sartorius.com)

### Poland

Sartorius Poland sp.z o.o.  
ul. Wrzesinska 70  
62-025 Kostrzyn

Phone +48.61.6473830  
Fax +48.61.6473839

### Russian Federation

LLC "Sartorius ICR" and LLC "Biohit"  
Uralskaya str. 4, Lit. B  
199155, Saint-Petersburg

Phone +7.812.327.5.327  
Fax +7.812.327.5.323

### Spain & Portugal

Sartorius Spain, S.A.  
Avda. de la Industria, 32  
Edificio PAYMA  
28108 Alcobendas (Madrid)

Phone Spain +34.902.123.367  
Phone Portugal +351.800.855.800  
Fax Spain +34.91.358.96.23  
Fax Portugal +351.800.855.799

### Switzerland

Sartorius Mechatronics Switzerland AG  
Ringstrasse 24a  
8317 Tagelswangen (ZH)

Phone +41.44.746.50.00  
Fax +41.44.746.50.50

### U.K.

Sartorius UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9QQ

Phone +44.1372.737159  
Fax +44.1372.726171

### Ukraine

LLC "Biohit"  
Post Box 440 "B"  
01001 Kiev, Ukraine

Phone +380.44.411.4918  
Fax +380.50.623.3162

## Americas

### USA

Sartorius Corporation  
5 Orville Drive, Suite 200  
Bohemia, NY 11716

Phone +1.631.254.4249  
Toll-free +1.800.635.2906  
Fax +1.631.254.4253

### Argentina

Sartorius Argentina S.A.  
Int. A. Avalos 4251  
B1605ECS Munro  
Buenos Aires

Phone +54.11.4721.0505  
Fax +54.11.4762.2333

### Brazil

Sartorius do Brasil Ltda  
Avenida Senador Vergueiro 2962  
São Bernardo do Campo  
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900  
Fax +55.11.4362.8901

### Canada

Sartorius Canada Inc.  
2179 Dunwin Drive #4  
Mississauga, ON L5L 1X2

Phone +1.905.569.7977  
Toll-Free +1.800.668.4234  
Fax +1.905.569.7021

### Mexico

Sartorius de México S.A. de C.V.  
Circuito Circunvalación Poniente  
No. 149  
Ciudad Satélite  
53100, Estado de México  
México

Phone +52.5555.62.1102  
Fax +52.5555.62.2942

## Asia | Pacific

### Australia

Sartorius Australia Pty. Ltd.  
Unit 5, 7-11 Rodeo Drive  
Dandenong South Vic 3175

Phone +61.3.8762.1800  
Fax +61.3.8762.1828

### China

Sartorius Scientific  
Instruments (Beijing) Co., Ltd.  
33 Yu An Road, Airport Industrial Park  
Zone B,  
Shunyi District, Beijing 101300,  
P.R.China

Phone +86.10.8042.6300  
Fax +86.10.8042.6486

### Hong Kong

Sartorius Hong Kong Ltd.  
Unit 1012, Lu Plaza  
2 Wing Yip Street  
Kwun Tong  
Kowloon, Hong Kong

Phone +852.2774.2678  
Fax +852.2766.3526

### India

Sartorius Weighing India Pvt. Ltd.  
#69/2-69/3, NH 48, Jakkasandra,  
Nelamangala Tq  
562 123 Bangalore, India

Phone +91.80.4350.5250  
Fax +91.80.4350.5253

### Japan

Sartorius Japan K.K.  
4th Fl., Daiwa Shinagawa North Bldg.  
8-11, Kita-Shinagawa 1-chome  
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.3740.5408  
Fax +81.3.3740.5406

### Malaysia

Sartorius Malaysia Sdn. Bhd  
Lot L3-E-3B, Enterprise 4  
Technology Park Malaysia  
Bukit Jalil  
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622  
Fax +60.3.8996.0755

### Singapore

Sartorius Singapore Pte. Ltd  
1 Science Park Road,  
The Capricorn, #05-08A,  
Singapore Science Park II  
Singapore 117528

Phone +65.6872.3966  
Fax +65.6778.2494

### South Korea

Sartorius Korea Ltd.  
8th Floor, Solid Space B/D,  
PanGyoYeok-Ro 220, Bundang-Gu  
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700  
Fax +82.31.622.5799

### Thailand

Sartorius (Thailand) Co. Ltd.  
129 Rama 9 Road,  
Huaykwang  
Bangkok 10310

Phone +66.2643.8361-6  
Fax +66.2643.8367



◀ [www.sartorius.com](http://www.sartorius.com)