

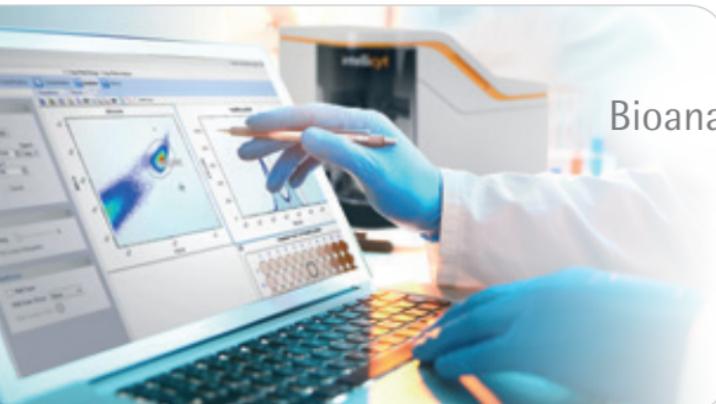


sartorius

Highlights of Our Lab Solutions

Issue Q4.2017 – Q1.2018





Bioanalytics



Bottle Closure &
Aseptic Fluid Transfer



Lab Filtration



Microbiological Analysis



Liquid Handling

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 Water



Lab Weighing



Centrifugation



Service & Training





iQue Screener PLUS

Get Cellular Insight at the Speed of Light



Perform rapid, multiplexed analysis of cells and beads in suspension with iQue Screener PLUS. 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

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



Derive deeper and more physiologically relevant information about your cells – plus real-time kinetic data with IncuCyte® S3.

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





MYCAP®

Make Aseptic Fluid Transfer in Bottles Possible



Robust, one-piece bottle closure with integral tubing for aseptic fluid transfer in bottles. MYCAP® is available on most bottles, flasks and centrifuge tubes. Easily and cost-effectively configure MYCAP® to meet specific requirements.

Features

- One-for-all silicone closure regardless of container type or tubing configuration
- One-piece bottle cap with integral tubing for aseptic fluid transfer
- Validated microbial barrier

Benefits

- Control quality and reduce validation costs
- Highly configurable in order to match MYCAP® to process requirements
- No risk of batch-to-batch or product-to-product contamination

Directly Harvest Your Mammalian Cell Cultures



The Sartoclear Dynamics® Lab P15 kit combines a syringe pre-filled with a filter aid for clarification and a 0.2 µm polyethersulfone filter for sterile filtration of up to 15 mL mammalian cell cultures all in a single step.

Features

- Cell culture filtration of up to 15 mL
- Sterile syringe filter with a prefilter and a 0.2 µm PES membrane
- Syringe pre-filled with 0.5 g of diatomaceous earth; includes a filling tube

Benefits

- No centrifugation step is needed
- Saves time due to higher flow rates
- No clogging of the membrane – one sterile filter per filtration run



Solvent Preparation



Arium® mini plus UV

- Up to 10 liter of constantly ultrapure water on demand
- Constantly low TOC content
- Constantly low conductance
- Bagtank technology to avoid contaminations

Creation of Standards

Cubis® individual with Q-App

- Fully automated preparation of reference standards
- CFR21 Part 11 compliant records
- Automatic solvent volume adjustment according to sample weight



Liquid Handling

Tacta® Mechanical Pipettes, Low-Retention Tips and Bottle

Top Dispensers

- Easy to clean
- High chemical resistance
- Optimum recovery of precious samples
- Reliable volume setting



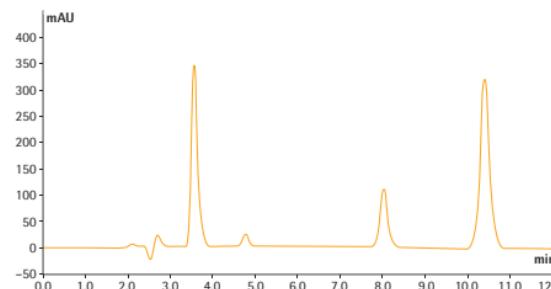
Filtration

Claristep® System or Minisart®

- Easy manual operation
- Filtrate up to 8 samples simultaneously with Claristep®
- Reliable removal of particles
- Syringe filters for different sample types



Clean HPLC Results



- Fewer ghost peaks
- Less background noise
- No column clogging
- Reliable results

HPLC – Clean Samples = Clean Results

HPLC as one of the most common analysis methods always pursues the target to gain reproducible and specific results. The quality of HPLC results is directly correlated to the quality of sample preparation. As this is demanding and time-consuming we at Sartorius support you to optimize your workflows with our solutions. By offering products for different HPLC sample preparation steps you can always rely on your HPLC results.





Microsart® Bacteria

Better Be on the Safe Side Fast!



Microsart® Bacteria detects >95 % of all known bacteria within a cell culture in only 3 hours. It is the perfect solution for rapid bacterial contamination control within regenerative medicine and bio-pharmaceutical in-process testing.

Features

- Internal control DNA is included in ready-to-use master mix
- Highly specific TaqMan® probes
- Non-infectious validation standards

Benefits

- The revolution in regenerative medicine – ensure ATMPs are bacteria-free prior to treatment
- No more worries about false positives – instead, obtain reliable results within 3 hours
- Rapid bioburden control during your whole biopharmaceutical process

Dispensing Made Easy



Prospenser & Prospenser Plus are safe, convenient and customizable. This makes them ideal for a wide range of dispensing needs.

Features

- The ceramic piston and all parts in the liquid pathway are highly resistant to chemicals
- The incremental volume setup ensures accurate and repeatable results
- Wide range of accessories available

Benefits

- Compatible with wide range of liquids
- Fully autoclavable for easy decontamination
- Easy and exact volume setting every time





arium[®] mini

Tailored to Your Specific Requirements



arium[®] mini lab water system best suited for your sensitive analytical application.

Features

- Innovative bagtank technology
- Multi-color touch-activated display with intuitive menu navigation
- Minimal footprint; only 28 cm wide

Benefits

- Closed system prevents secondary contamination
- Easy to operate without a user manual, even when you wear gloves
- Saves valuable space anywhere in your lab

Forget the Inconvenience of Transferring Samples



Avoid transfer errors and sample loss: reliably weigh small samples directly into heavy sample containers of up to 50 mL with Cubis® High-Capacity Micro Balances.

Features

- Flexible and adjustable sample holders
- The world's highest resolution in a micro weighing instrument with up to 60 million divisions
- Weighing capacity of more than 60 g and a readability of 1 µg

Benefits

- Ergonomic and reliable dosing
- Simplified process because no sample transfer is necessary
- No more wasting of expensive samples because the Cubis® mirco balance enables a low sample weight





Centrisart® G-26C

Centrifuge at
High Speed
While Refrigerating
Your Samples



The refrigerated centrifuge Centrisart® G-26C achieves a gravitational field of up to 62,000 g at a maximum speed of 26,000 rpm.

Features

- Easy to operate; patented one-knob control
- Large LCD display
- Process library with program preview

Benefits

- Easy to operate thanks to intuitive menu
- All parameters on one easy-to-read screen
- No hassle with trying to find the right program

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



Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Germany



◆ www.sartorius.com