



High-Throughput
and Essential Lab Tools
to Streamline Cell Line
Development

Simplifying Progress

SARTORIUS

Innovative Technologies to Select the Right Clone



Gene Cloning and Initial Clone Selection
Rapid screening of minipools for product titer, cell health and growth



Clone Selection and Confirmatory Analysis
Early identification of high productivity clones with optimal critical quality attributes



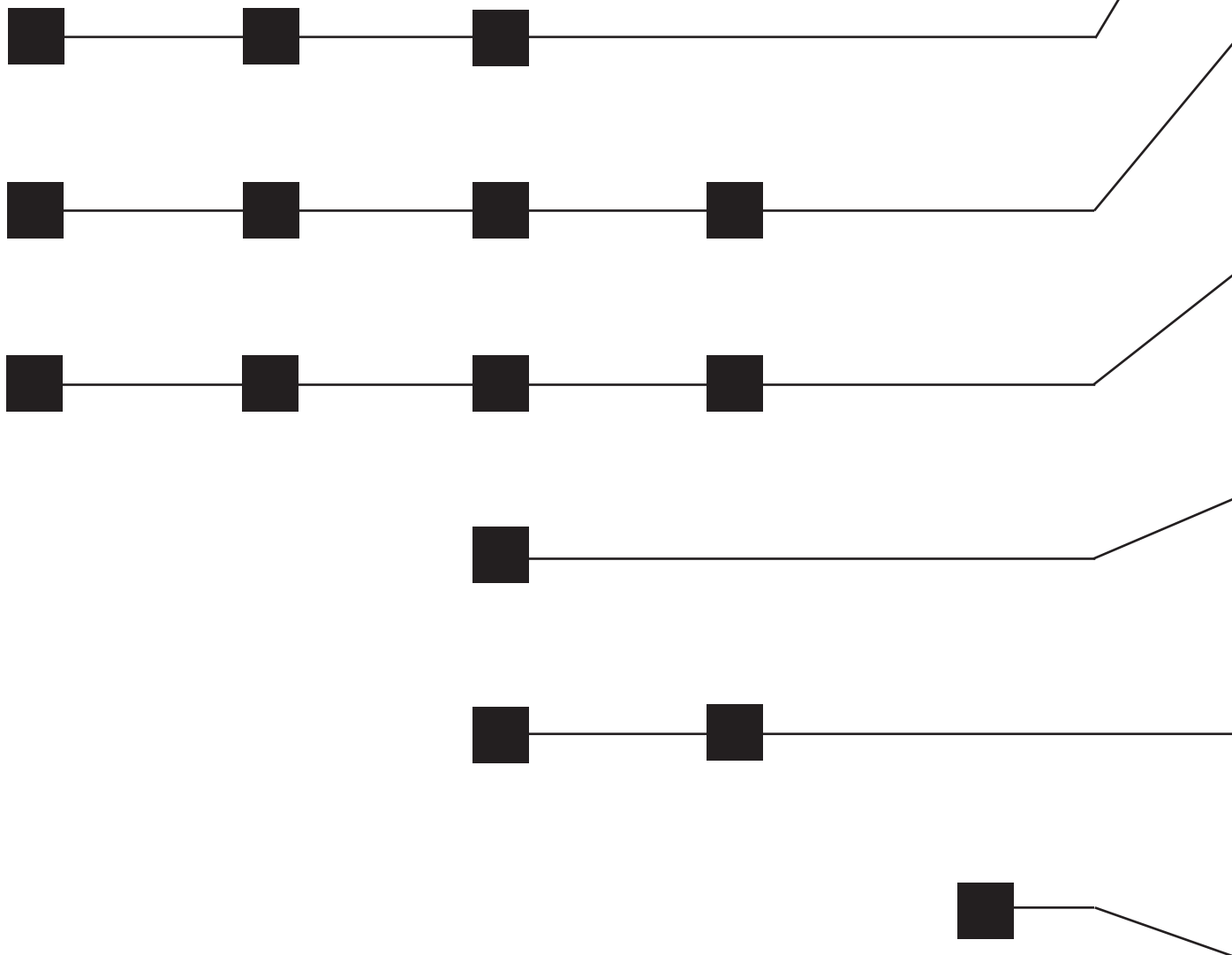
Cultivation and Media Optimization
Screen culture conditions and media formulations for optimal growth and protein production



Cell Line Evaluation and Characterization
Characterize top clone for stability, productivity, and quality attributes of the protein



Cell Banking
Research Cell Bank (RCB), Master Cell Bank (MCB), and Working Cell Bank (WCB) preparation



Solutions & Services for Cell Line Development

Select and screen single cells for antibody discovery

The CellCelector single cell and colony picking platform is used for high-throughput single cell cloning, allowing fast generation of clonal production cell lines. Cell lines are produced in one round while providing in-process, image-verified monoclonality proof.

- Single cell cloning and picking
- Integrated monoclonality, viability and productivity proof
- Characterize tens of thousands of clones in parallel

CellCelector Flex Automated Cell Imaging and Picking System



Perform protein quantitation and kinetics characterization assays

Measure molecular interactions in real-time without the need for detection reagents and enable advanced, fast, and fluidics-free direct detection of specific biomolecules.

- Rapidly determine product titer
- Screen for specific glycosylation content
- Perform off-rate analysis using purified and unpurified samples

Octet® Bio-Layer Interferometry (BLI) Detection Systems for Label-Free Detection



Multiparametric and actionable results for more insight

A high-throughput, suspension cell and bead-based analysis platform for antibody screening and functional profiling. The fastest way to generate high content data from small sample volumes.

- Screen your libraries in days, not months
- Sample volumes as low as 1 µL, reduce reagent costs
- Screen cells in suspension and targets in their natural conformation

iQue® Advanced Flow Cytometry Platform



Improve the overall efficiency and speed of your cell line and medias

Ambr® 15 Cell Culture is a high-throughput automated micro bioreactor for predictive cell line and media screening. Rapidly evaluate multiple bio reactor cultures at microscale, thus increasing your productivity in cell line development and saving on materials and labor.

- Rapid screening of clones and media optimization
- Data-rich culture profiles with automated sampling
- Large multifactorial experimental designs (DoE) in real-time

Ambr® 15 Cell Culture Bioreactor System



An automated, high throughput benchtop system that performs a one-step clarification and purification process to streamline and automate mAb cell culture harvest coming from Ambr® 15 and other suspension cell cultures.

- Predefined templates, flexible process selection, and error recovery creates a walk-away process that reduces downtime
- Saves space and simplifies workflows by removing the need for separate centrifugation, filtration, and chromatography systems
- Detailed data recording and auditing for extended data analysis capabilities

StreamLink® CC 15 mAb Sample Preparation



Produce large, homogenous banks of cells under cGMP conditions

Fill-it Automated Cell Banking System offers improved consistency and shortens process times compared to manual processing, giving the potential to increase batch sizes for cell banking and dramatically reduce QC costs.

- Fits in a standard biosafety cabinet
- Facilitates easy loading and operation
- De-caps racks of cryovials held in either 96-, 48- or 24-vial racks
- Provides automated cell banking compatible with wide microtube range

Fill-it Automated Cell Banking System



Lab Essentials for Cell Line Development

Everyday lab essential products play a critical foundational role in the success of cell line development and are invaluable to lab technicians and scientists. Basic tasks like pipetting, filtration and purification of small volumes can be time-consuming and tedious. Sartorius' lab essentials help streamline routine tasks and achieve more accurate results by:



Increasing reproducibility of your results



Avoiding contamination in your processes



Increasing safety and reduce fatigue with ergonomic and easy-to-use products

Pipetting

Picus® 2 electronic pipettes ensure accurate results with an unbeatable ergonomic design. Connect Picus® 2 to your mobile device to smoothly run sample preparation workflows and adjust settings automatically, taking your productivity to the next level.



Midi Plus are lightweight, easy-to-use pipette controllers for large volumes applications.



Standard and Buffer Preparation

The **Cubis® Dosing System** offers semi-automatic preparation and documentation of highly reproducible calibration standards for chromatography and spectroscopy.



Vivapure® and Sartobind® Membrane Adsorbers
Rapid, easy target purification and contaminant removal by anion or cation exchange chromatography.



Cubis® II Laboratory Balances

100% Configurable Laboratory Balance

With completely configurable hardware, software and connectivity, Cubis® II offers a high-performance balance that will align with your unique demands and compliance requirements.



Arium® Type 1 Ultrapure Water Systems

ASTM Type 1 water, for general, analytic and life science applications, ensuring consistent, high-quality results, at all times.



Mycoplasma Detection

Keep your cell cultures contaminant-free and compliant with **Microsart® Mycoplasma qPCR kits**.



Lab Instrument Services

Certified and accredited services. Including:

- Installation/Equipment qualification (IQ/OQ)
- Extended warranty
- Calibration
- Preventive maintenance
- Service contracts, original spare parts



Clarification, Concentration and Purification of Proteins

Sartoclear Dynamics® Lab



Vivaspin® and Vivaflow® Ultrafiltration Devices



Vivapure® and Sartobind® Membrane Adsorbers



Minisart® Syringe filters Sartolab® RF | BT Vacuum Filtration Units



Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

UK

Sartorius Stedim Lab Ltd.
Sperry Way, Stonehouse Park
GL10 3UT Stonehouse,
Gloucestershire, UK
Phone +44 1453 821972

 For further contacts, visit:

www.sartorius.com