



# Bioprocessing Media and Buffers

Solutions That Adapt to  
Your Needs

Simplifying Progress

**SARTORIUS**





## Optimization of Manufacturing Processes

### Maximizing Efficiency & Reducing Time-to-Market

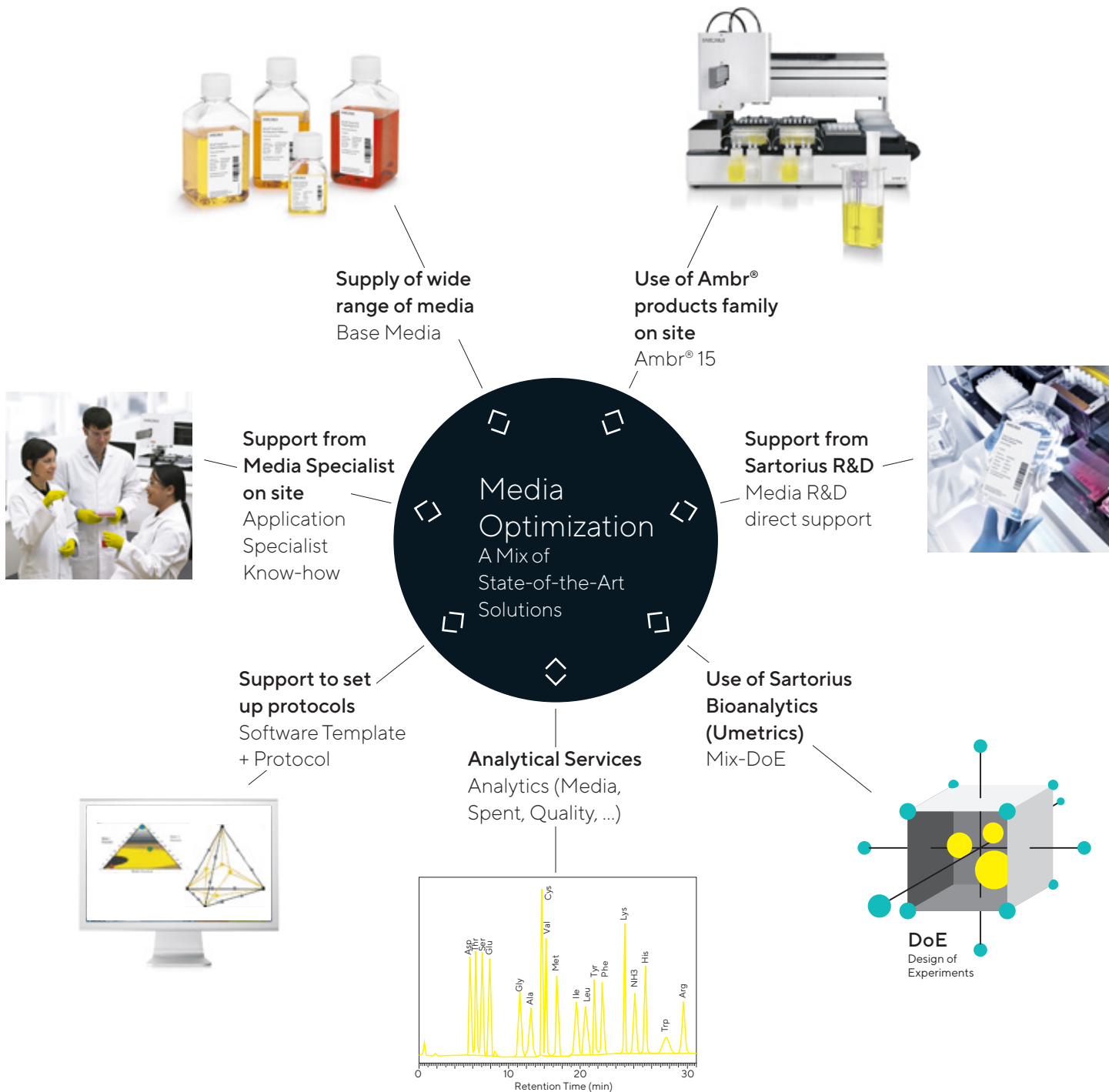
With the growing needs of globally accessible therapeutics and competitive pressure in the biopharmaceutical market, the industry is challenged to find innovative treatments for a wider selection of indicators.

With the number of parameters impacting the manufacturing process and its output, supported systems approaches are more and more becoming an integrated necessity in process development.

The media and buffers offered by Sartorius are an essential component in this optimization with a significant role in shaping the overall manufacturing process.

### Benefit From Our Media Expertise & Media Optimization Service

- Experienced in Process Optimization and Buffer usage – Only a few steps from you.
- Wide range of media formulation commercially available
- Design Media and Feeds suitable for your needs



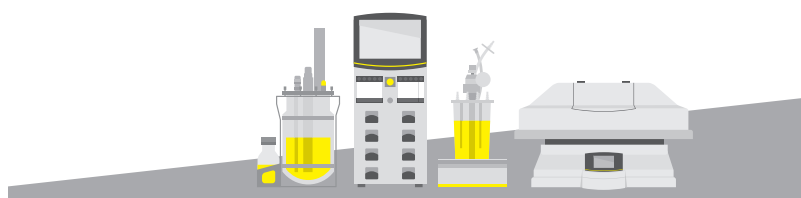


## Optimizing Your Processes and Media Composition

Taking into account the process and all variable elements included will significantly improve the development stages and time-to-market. Process parameters, cells and cell culture media interact in a challenging way to define the productivity and quality of the protein output. Process-integrated media service provided by Sartorius will demonstrate the benefits it can bring to you.

## Scalability

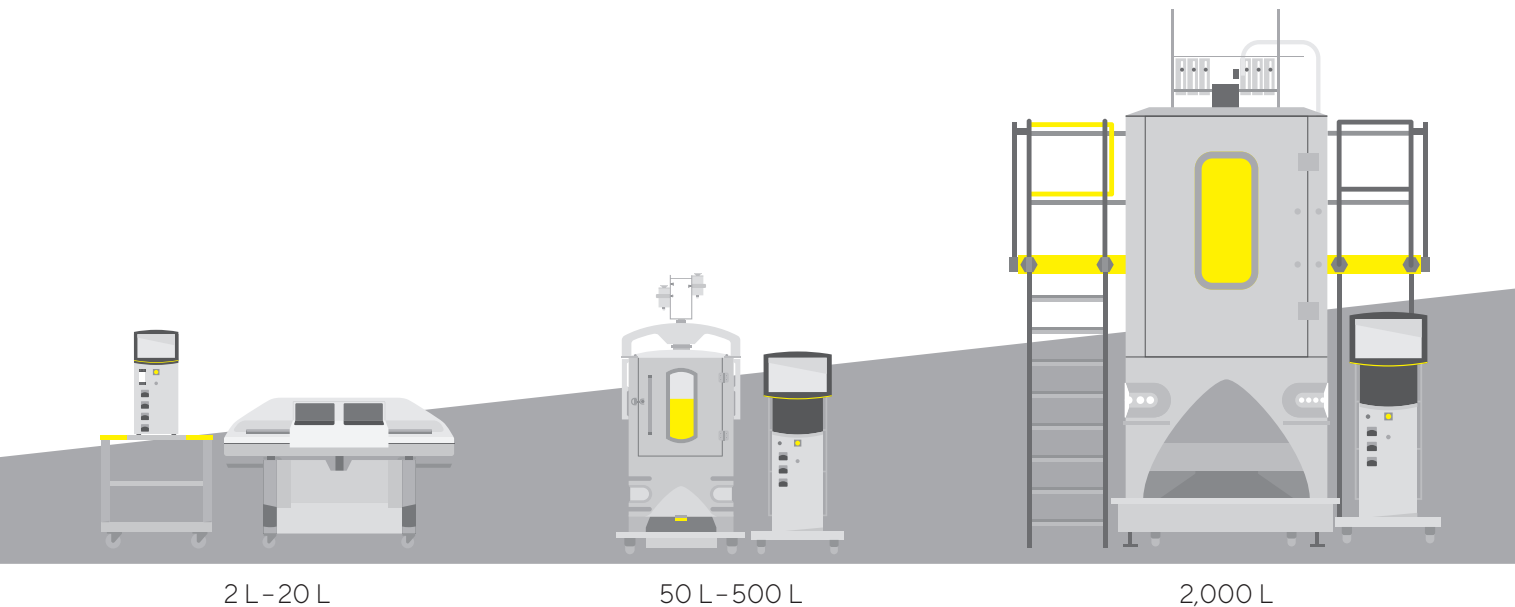
Scalability from early process development to commercial production



250 mL – 10 L

Your Goals	Your Needs	Our Solution
Faster and smart way to develop your process	Quick Support while improving your processes & Benefits of good and structured media optimization	Set up your process parameters quicker with dedicated expert know-how and availability with the help of smart tools.
Better knowledge of your process needs and parameters. Getting optimized media and feeds according to your requirements.	Improvement of your media composition & cell-based process	Our Media Specialists help you analyze data and set up the optimal recipe focusing on accelerating development time and transfer to production.
Getting Media and Buffer Products in the best quality and suitable quantity according to the step in the manufacturing process.	Wide Choice of formats and primary packaging for pilot scale Cell Culture Media and Buffers and further manufacturing use in GMP	Committed Project Management dedicated to handle your inquiry from RFQ to delivery. Manufacturing sites capable to deliver products in liquid and powder in different packaging formats





### Your Goals

Testing new formulations within process development stage. Getting small volumes fast with same raw material quality than at large scale (suitable for GMP usage)

Getting Media and Buffers Products in optimal packaging for your needs. Optimising your manufacturing COGs.

Business continuity

### Your Needs

Pilot Scale Cell Culture Media and Buffers: Fast & Clean

Production of Media & Buffers for further manufacturing use in GMP in wide choice of formats and primary packaging

Solid partner

### Our Solution

Dedicated Project Manager to handle your custom inquiry from RFQ to delivery. Two Manufacturing sites

Two manufacturing sites capable to deliver products in liquid and powder in different packaging formats

Sartorius: Plural Manufacturers with long-term agreements



# Delivering Your Media & Buffers

## High Quality Manufacturing Capabilities

- ISO-Certified manufacturing sites on different continents
- Liquid and Powder production
- Capable of transferring production from one site to another to increase risk mitigation
- Same high-quality of raw material used for pilot production | batch production

Pilot Capabilities		Large Capabilities per Batch (according to cGMP)	
Liquid	Powder	Liquid	Powder
1-100 L	1-20 kg	Up to 10,000 L	Up to 7,000 kg

## Ensured Operational Excellence

- Media manufacturing facilities have multiple environmental monitoring and cleaning protocols in place to ensure operational excellence.
- ISO Certified
  - Site wide environmental monitoring
  - Incoming Goods control
  - Wide Variety of QC tests for Release
  - Documentation for Regulatory Dossier



Powder Media | Buffers for Large Scale Manufacturing Use

Liquid Media or Buffers for Small to Large Scale Manufacturing Use

Safe & stable secondary packaging solution for Media or Buffers



## Full Range Media – Your Complete Solution

### Off-the-Shelf or Custom

- Wide range of Cell Culture Media and Buffers
- Media Optimization (Support for screening of commercially available standard offerings or personalized dedicated support)
- Fast and Clean Pilot Offering (Expedited production option for liquid and powders using same raw material sourcing as our large scale manufacturing)

### Powder Media

- Powder media manufacturing techniques for optimal product (High solubility, batch consistency and performance)

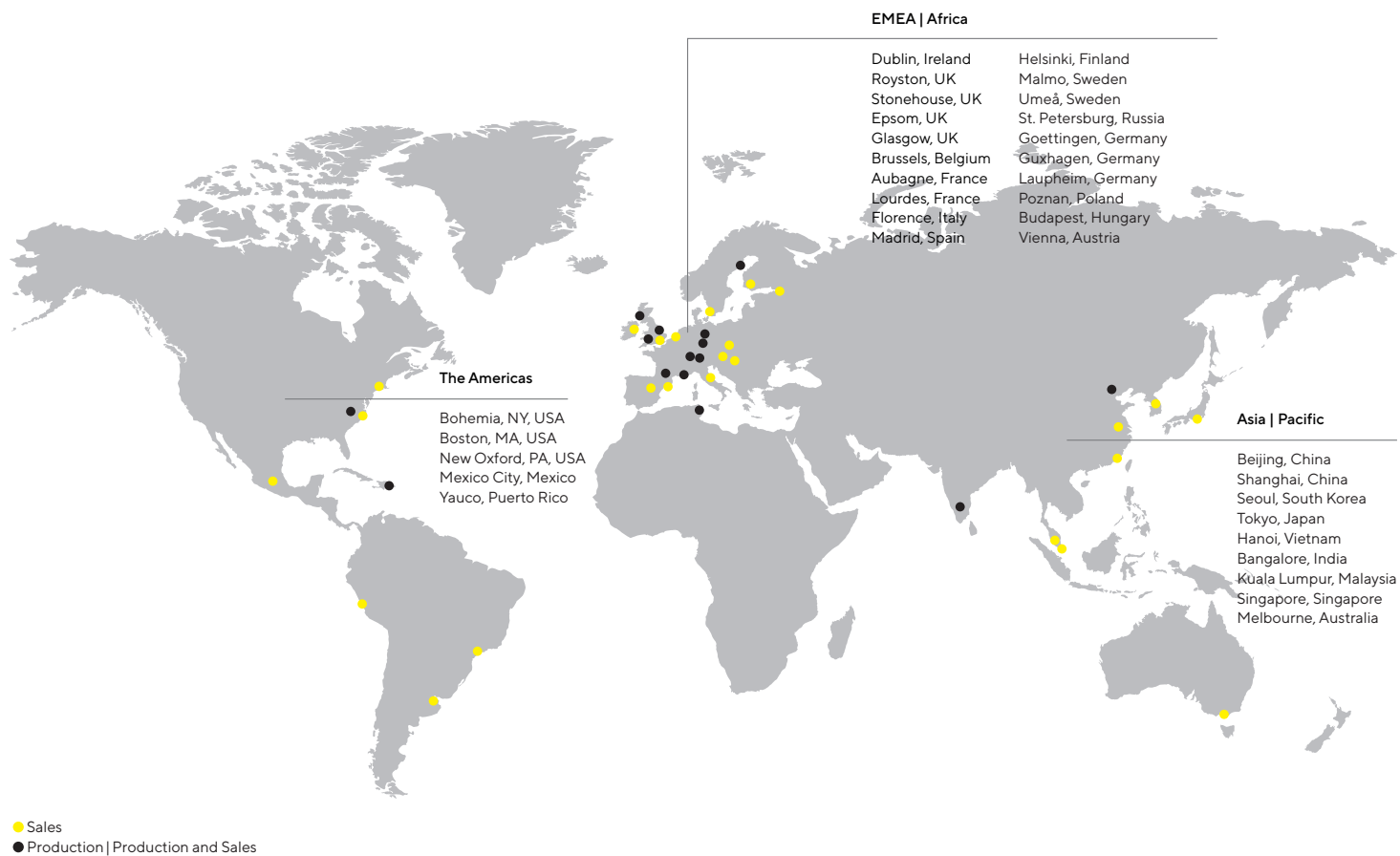
### Buffers

- Available in a range of sizes as sterile-filtered liquids or in modern pre-measured powder formats

## Everything at One Glance

Benefits   Your Need		Features   Our Solution
Media and buffers in Sartorius Stedim bags	>	Qualified Film and Assurance of Supply
All Media Filtered through Sartopore® 2 family of proprietary filters	>	High level of Scale Up success and quality
Pre-weighed Powder Media and Buffers in bags and liners	>	Innovative and cost-effective solution for media preparation
Products fully integrated into the Sartorius upstream and downstream solutions	>	One single point of contact for all project needs

# Sartorius Stedim Biotech Facilities



## Germany

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

## USA

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

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

Specifications subject to change without notice.

© 2021 Sartorius Stedim France S.A.S., Avenue de Jouques - Zone Industrielle Les Paluds - C.S. 71058, 13781 Aubagne Cedex, France

Status: 06 | 02 | 2021