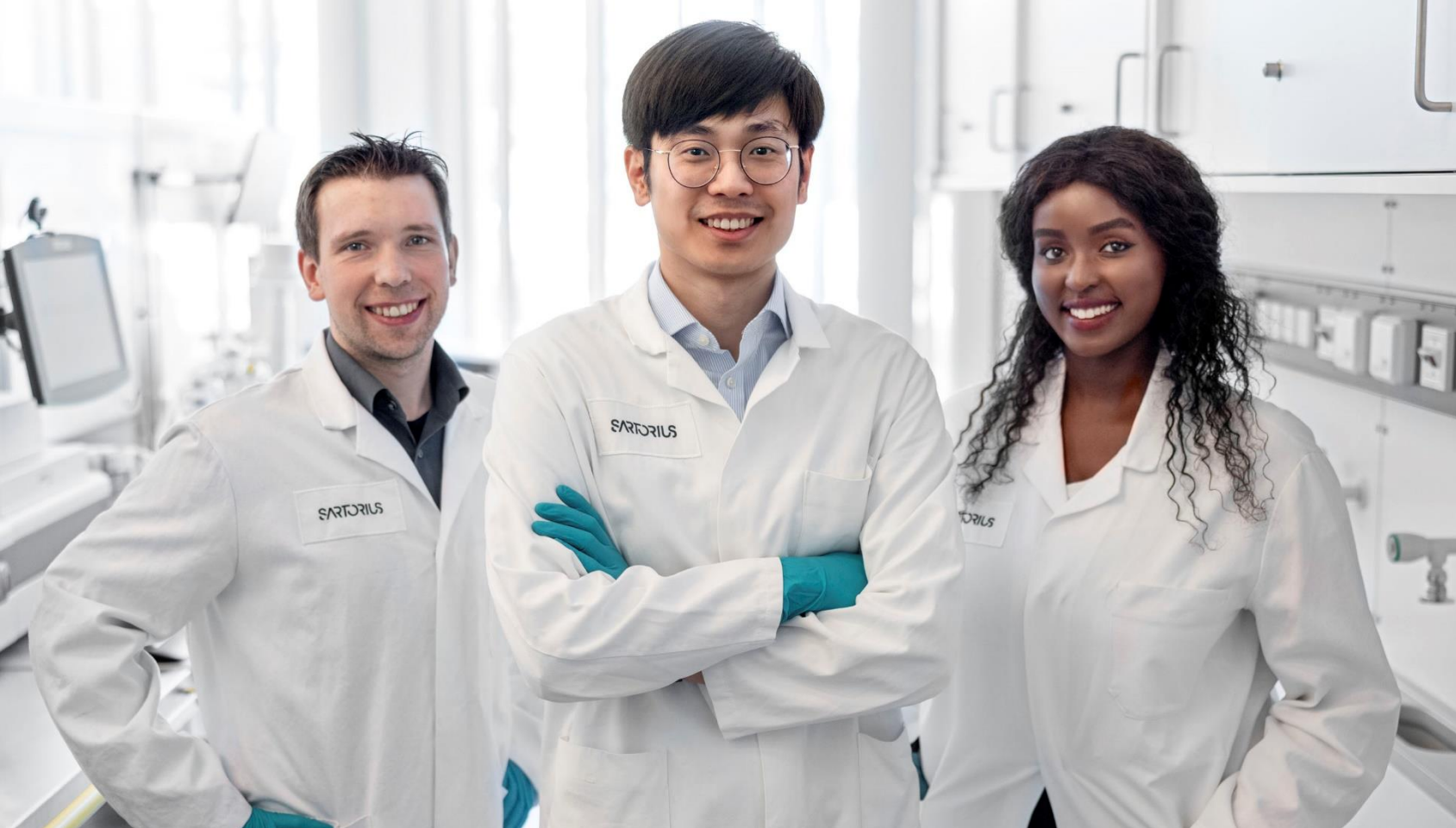




Simplifying Progress

Welcome to Sartorius

February 2025

SARTORIUS

Our ambition: Better health for more people

Mission

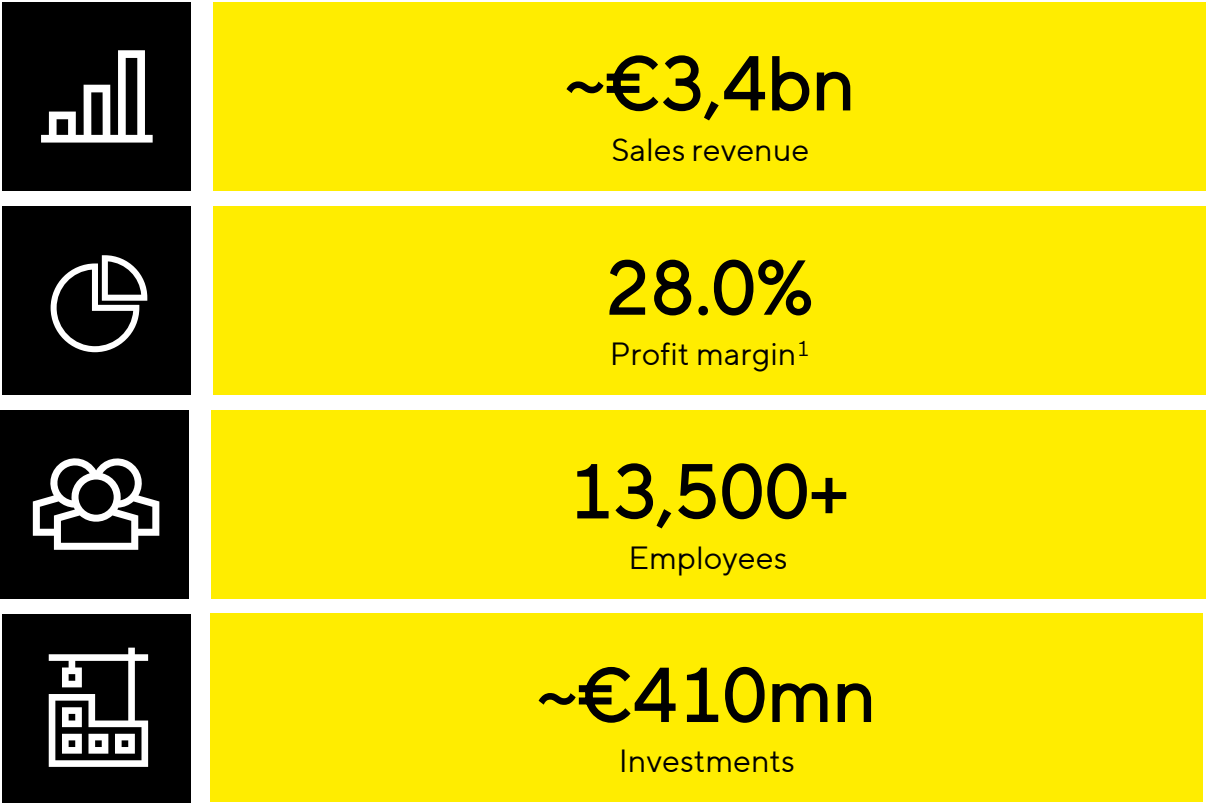
We empower scientists and engineers to simplify and accelerate progress in life science and bioprocessing, enabling the development of new and better therapies and more affordable medicine.

Vision

We are a magnet and dynamic platform for pioneers and leading experts in our field. We bring creative minds together for a common goal: technological breakthroughs that lead to better health for more people.



Sartorius in 2024 numbers



1 Underlying EBITDA margin



13,500 employees united by strong values

Sustainability

Growing profitably and acting responsibly towards all stakeholders



Openness

Driving change and progress internally and externally



Enjoyment

Working in an energetic and rewarding environment



112

Nationalities

52%

Younger than 40 years

39%

Women

50%

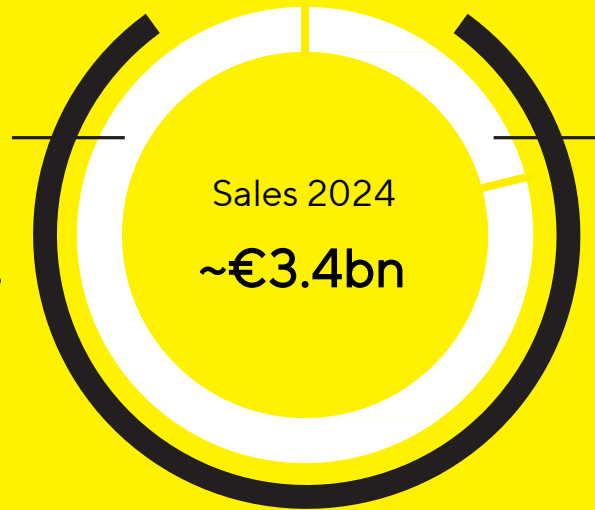
Joined since 2019

Two divisions with a clear focus on the life science industry

Offers a wide array of innovative technologies for the manufacture of biopharmaceuticals and vaccines as well as cell and gene therapeutics



80%
Bioprocess Solutions
Division



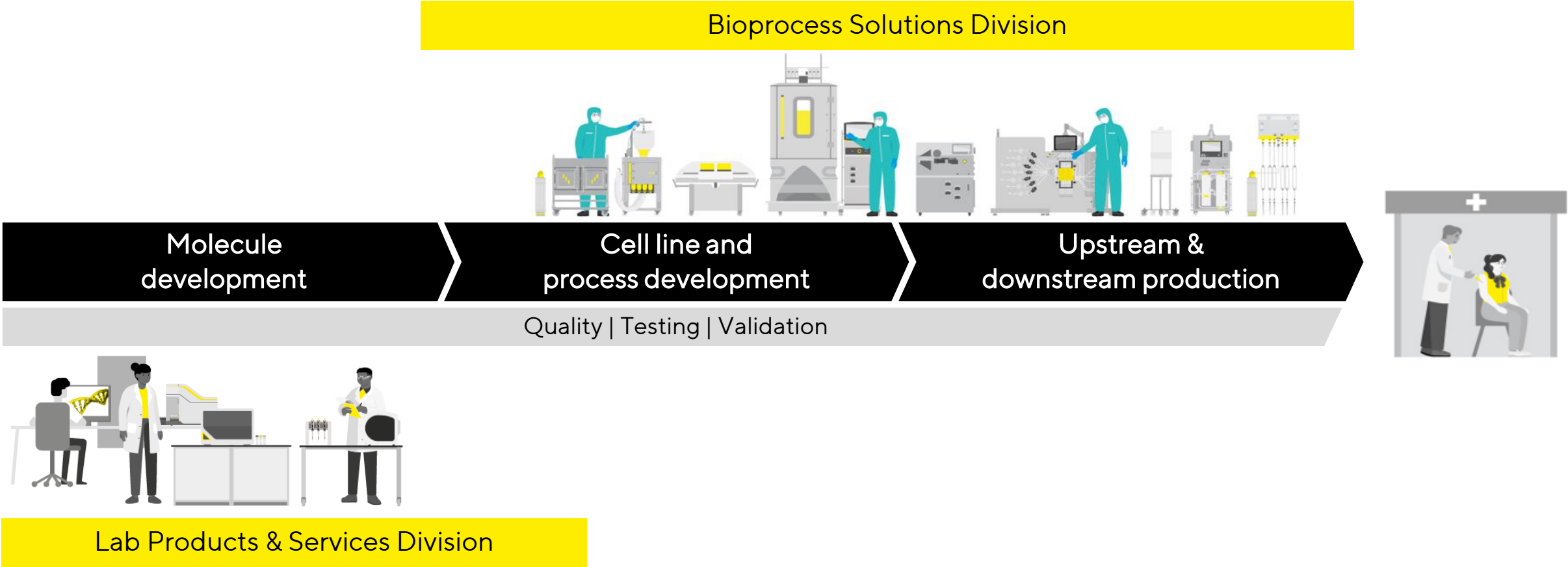
~85%
with life science
customers

20%
Lab Products & Services
Division



Specializes in
equipment and
technologies for life
science research and
pharmaceutical
laboratories

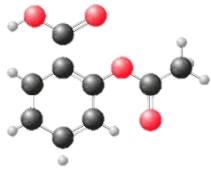
Simplifying Progress: Enabling the discovery and production of biopharmaceuticals



What are biopharmaceuticals?

Chemical drugs

Aspirin
($C_9H_8O_4$)



Small, simple
molecules



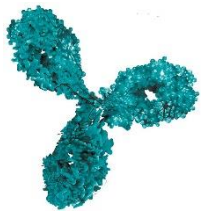
Obtained via
chemical synthesis



Administered
in various ways

Biopharmaceuticals

Monoclonal
antibody



Large molecules,
cells or viruses



Produced with living cells
in cell culture processes



Administered
mainly intravenously

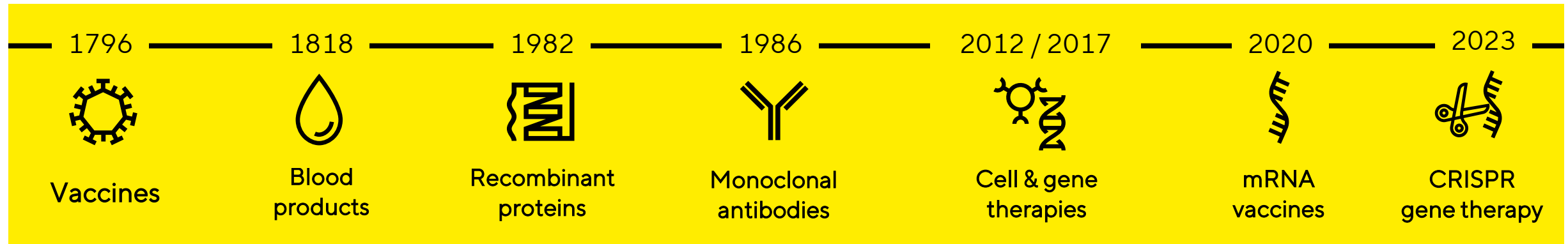
Advantages of biopharmaceuticals

- Target only diseased cells
- Fewer side effects
- Allow first-time or improved treatment of illnesses, e.g. cancer, multiple sclerosis, rheumatism



The promises of biologics: From established to new concepts

Breakthroughs at an ever-increasing pace and new forms of treatment offer hope for patients



19mn

Cancer diagnoses
per year

400mn

People with rare diseases,
>90 % without treatment options

1 in 9

People >65 years suffer
from Alzheimer's disease

Many biotech medications are highly effective, but very expensive

Rheumatism

Monoclonal antibodies



~€12,000

Annual cost of treatment in GER

Blood cancer

CAR-T cell therapy



~€240,000

Cost per treatment in Europe

Leukodystrophy

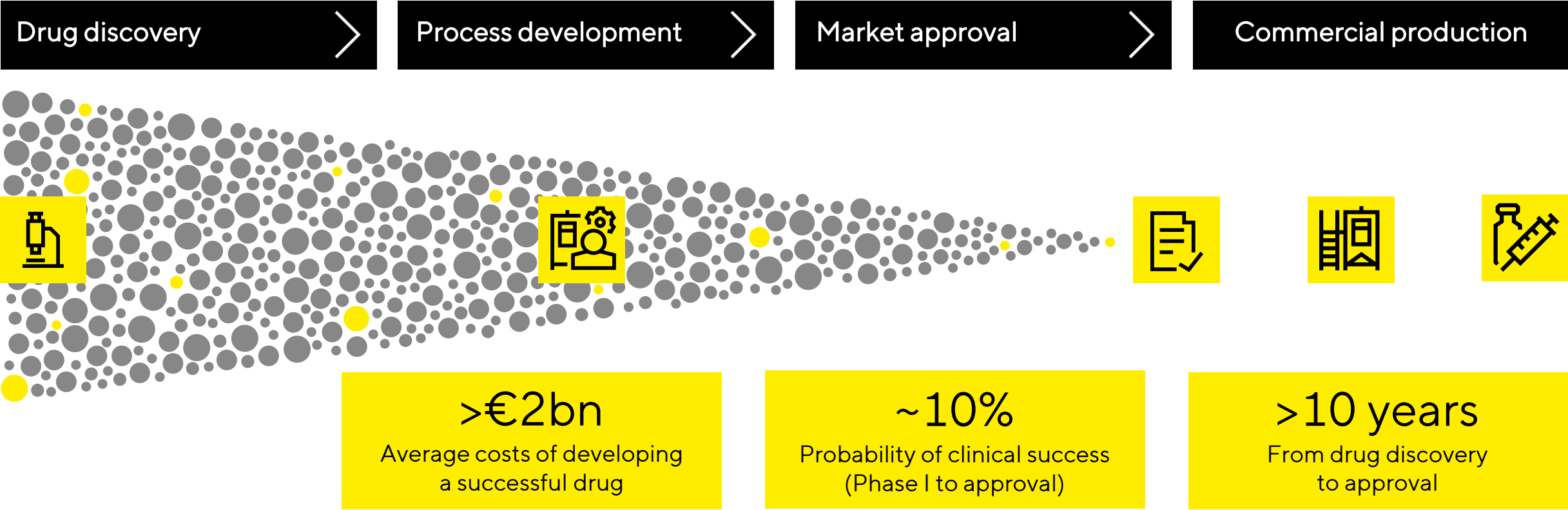
Gene therapy



~\$4.25mn

Cost per treatment in the USA

Only one out of 10,000 new drug candidates reaches the market



Schematic example of biologic drug discovery with data from the Association of the British Pharmaceutical Industry

Helping researchers in the labs to accelerate drug discovery



Our ambition
Reduce costly trial & error



Our solutions
Powerful bioanalytical tools and lab essentials



- ☒ Understand diseases
- ☒ Conduct experiments
- ☒ Make better use of data
- ☒ Identify molecules
- ☒ Develop new drugs



Live cell analysis



Cell characterization



Protein analysis



Cell selection



Pipetting



Weighing



Filtration



Water purification

Once discovered also the manufacture of biologics remains complex



Providing technologies to increase productivity in bioprocessing



Our ambition
More efficient bioprocessing



Our solutions
Technologies for all phases of drug manufacturing



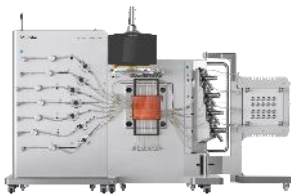
- ✓ Set up safe and flexible processes
- ✓ Increase resource and material efficiency
- ✓ Reduce investments
- ✓ Enhance product yield



Media & reagents



Bioreactors & fermenters



Chromatography



Filtration



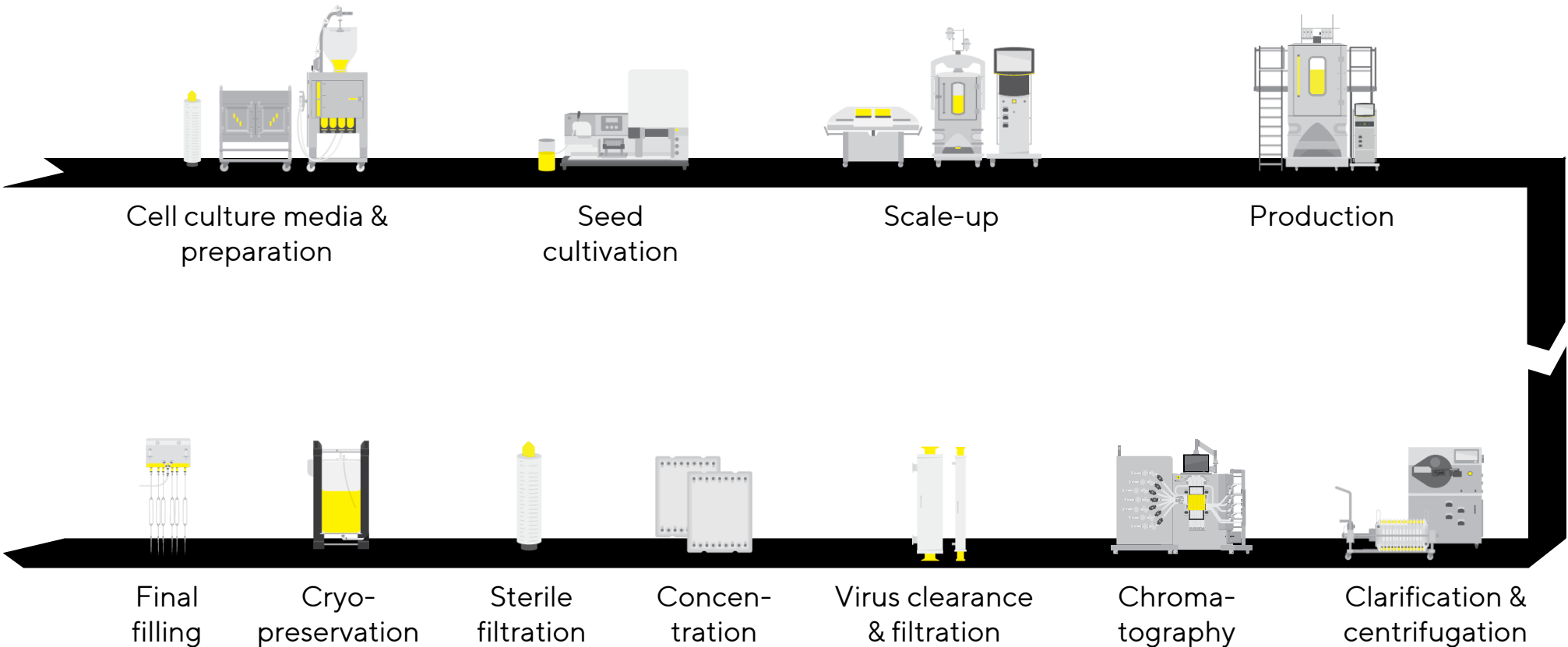
Fluid management



Data analysis

Offering integrated end-to-end manufacturing solutions

Upstream
Production of
the desired
drug



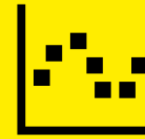
Helping customers to reduce their environmental footprint



Single-use
technologies



Process
intensification



Data
analytics



- No water-/energy-intensive cleaning
- Less cleanroom space, less energy usage
- Higher yields



Biostat® STR Bioreactors



Linkit® AX



Resolute® BioSMB

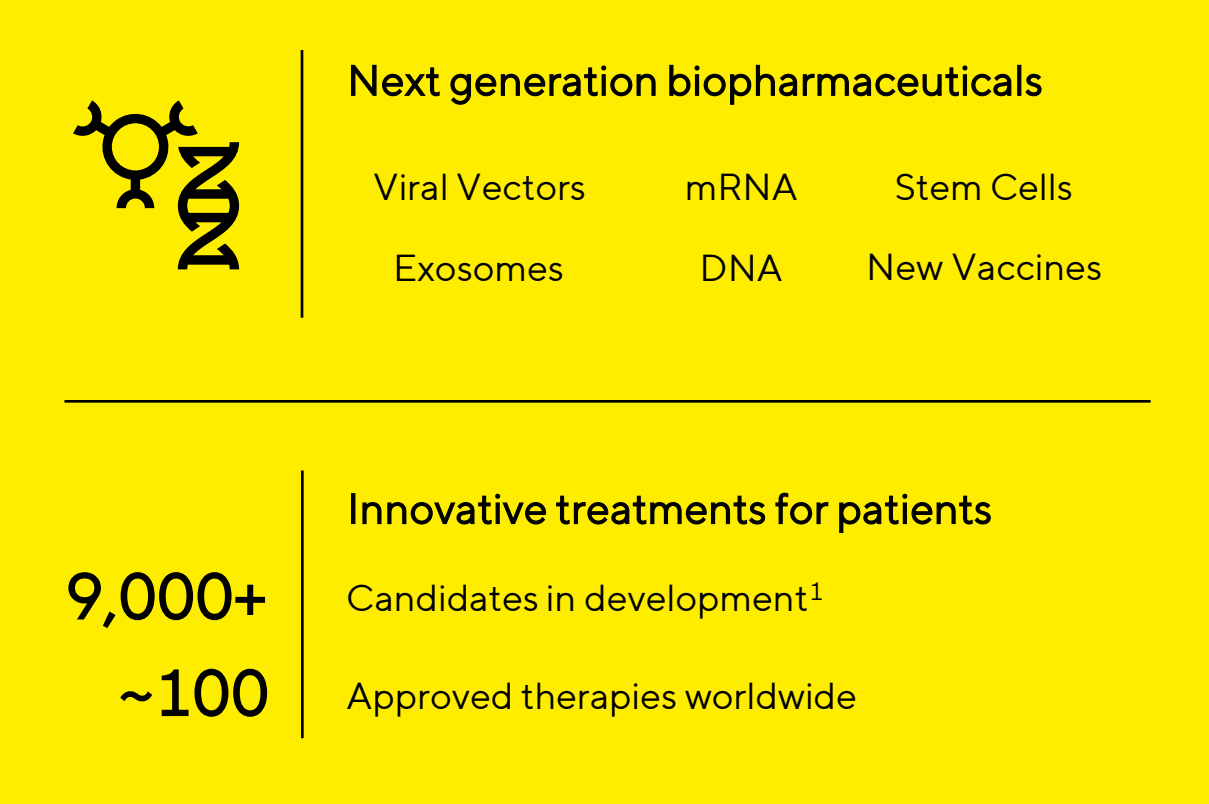


Ambr® 250HT Perfusion



Umetrics Suite

Cell and gene therapies: New modalities, new challenges

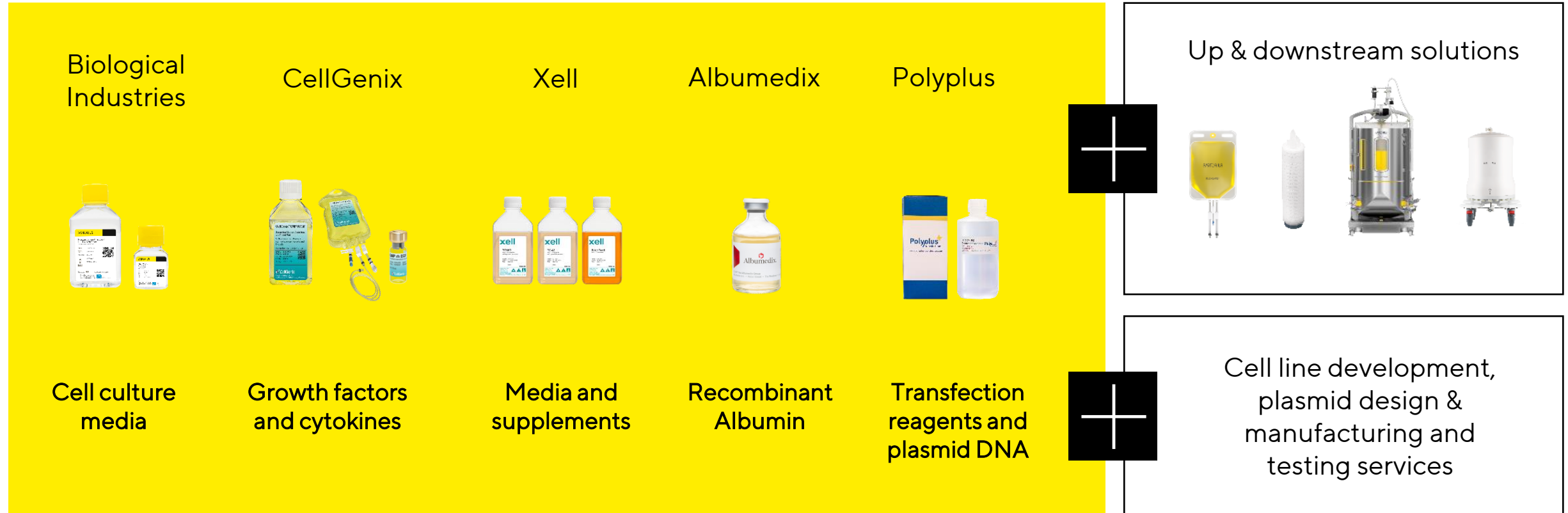


¹ Global Data, October 2024, CGT = Cell & Gene Therapies

This section features a black arrow-shaped banner at the top with a gear and upward arrow icon, followed by the text 'Enabling CGT: High need for innovation'. Below this, a list of five challenges is presented, each preceded by a plus sign. To the right of the list is an illustration of a scientist in a white lab coat holding a test tube.

- + Robustness of production process
- + Yields, productivity, cost-per-dose
- + Time-to-market
- + Regulation under development
- + Manufacturing close to patients

Extensive technology platform for cell and gene therapy applications



How we innovate



Own Product Development

in areas of core competencies

Separation

Cell cultivation
technologies

Lab analysis

Automation

Recent focus; selected projects



Acquisitions

of complementary technologies

Media & reagents

Chromatography

Bioanalytical instruments



Cooperations

with partners leading in their fields

Process
intensification

Bioprinting

Artificial
intelligence

Cell models

Resilient global setup, well prepared for further growth



Customer proximity



Global resilience



Redundant capacities



Operational excellence

~€410mn
Investments 2024

Expansion of production & R&D capacities; selected current projects

Opened in 2024

Selected projects



Marlborough, USA
Center for Bioprocess Innovation



Ann Arbor, USA
Bioanalytical products & services

Current investments

Selected projects



Göttingen, Germany
Filters



Songdo, South Korea
Media, bags, filtration, labs, training

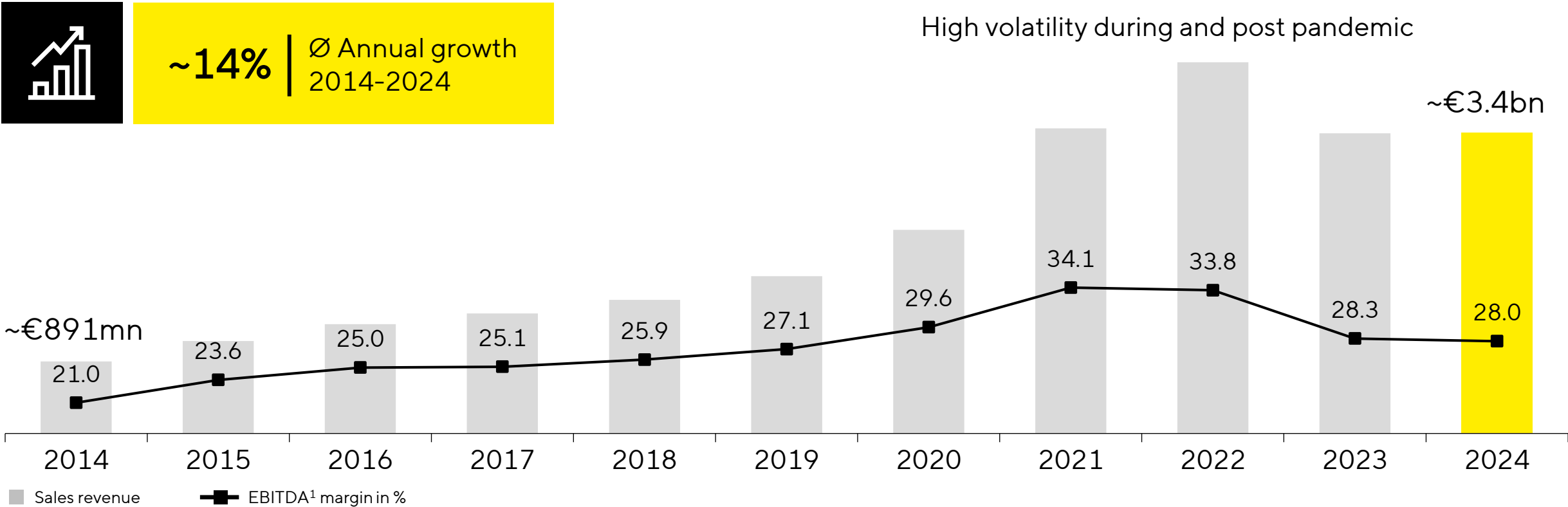


Freiburg, Germany
Cell culture components



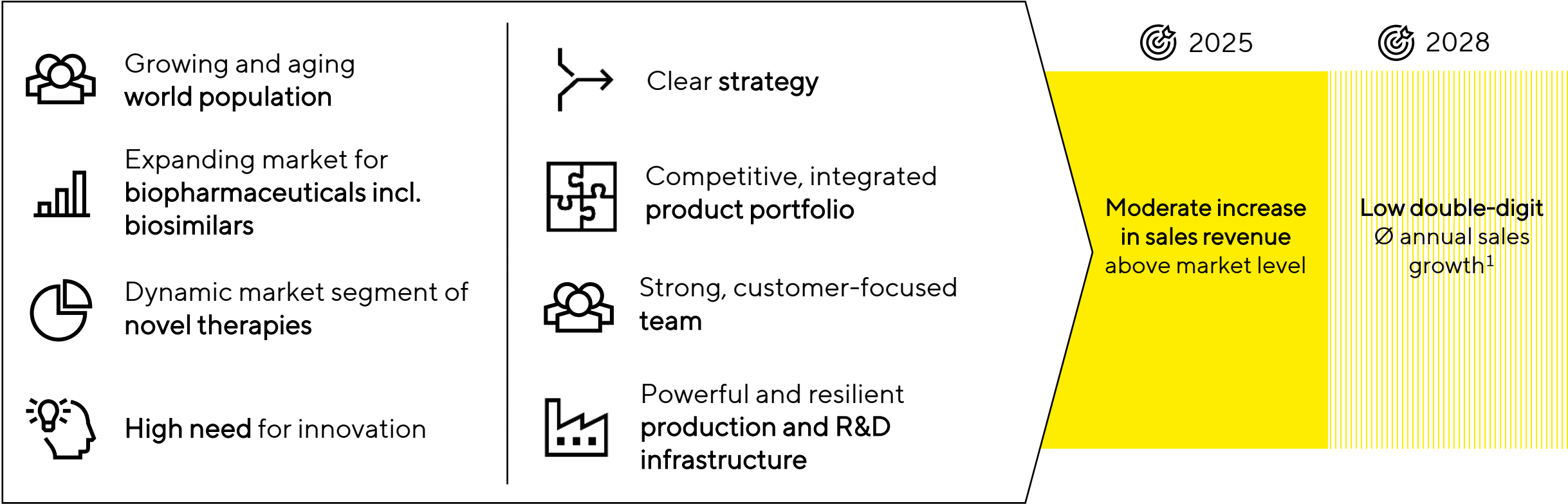
Aubagne, France
Clean rooms, product development, labs

Strong financial track record; profitable growth



Sales CAGR for continued operations, in constant currencies; EBITDA margin excluding extraordinary items

Ambitions backed by strong market fundamentals



1 Thereof around a fifth through acquisitions

Ambitious sustainability agenda in implementation

2030 mid-term ambitions

~10%

av. reduction of CO₂eq emission intensity p.a. (Scopes 1-3)

Zero

avoidable¹ emissions in Scopes 1 and 2

100%

electricity from renewable sources



¹ Process emissions from membrane manufacturing are currently considered unavoidable based on available technology

2045 long-term ambition

Net-zero emissions

Decarbonization of the entire supply chain in collaboration with suppliers and customers



Simplifying Progress

Sartorius AG
Corporate Communications & IR
Otto-Brenner-Straße 20
37079 Göttingen
Germany

SARTORIUS