

SARTORIUS

Simplifying Progress

Transformation of upstream bioprocessing in the digital era

World Bioprocessing Pharma 4.0 Summit- October 2021



Sartorius in brief

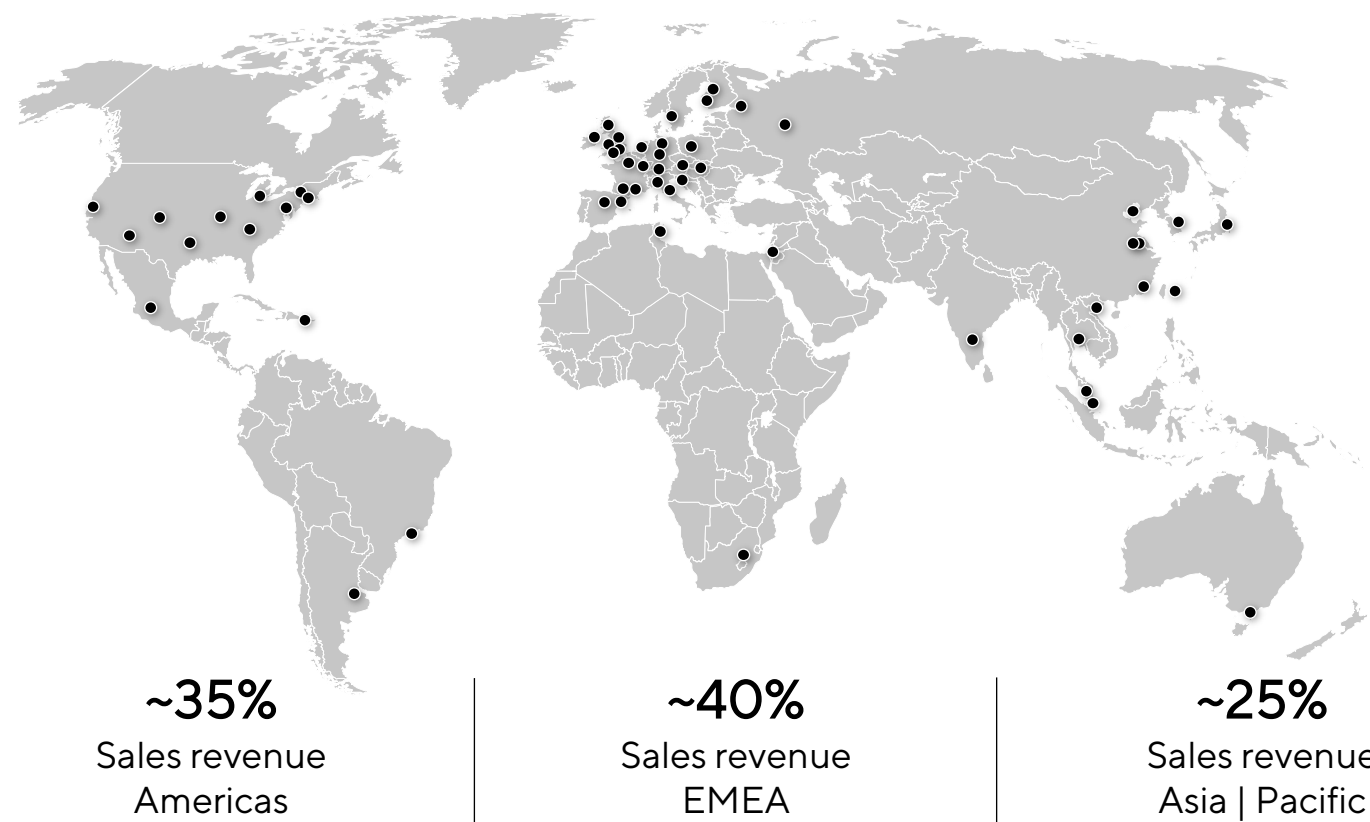
 **60+**
Locations worldwide,
headquartered in Göttingen, Germany

 **10,600+**
Employees¹

 **~€2.3bn**
Sales revenue¹

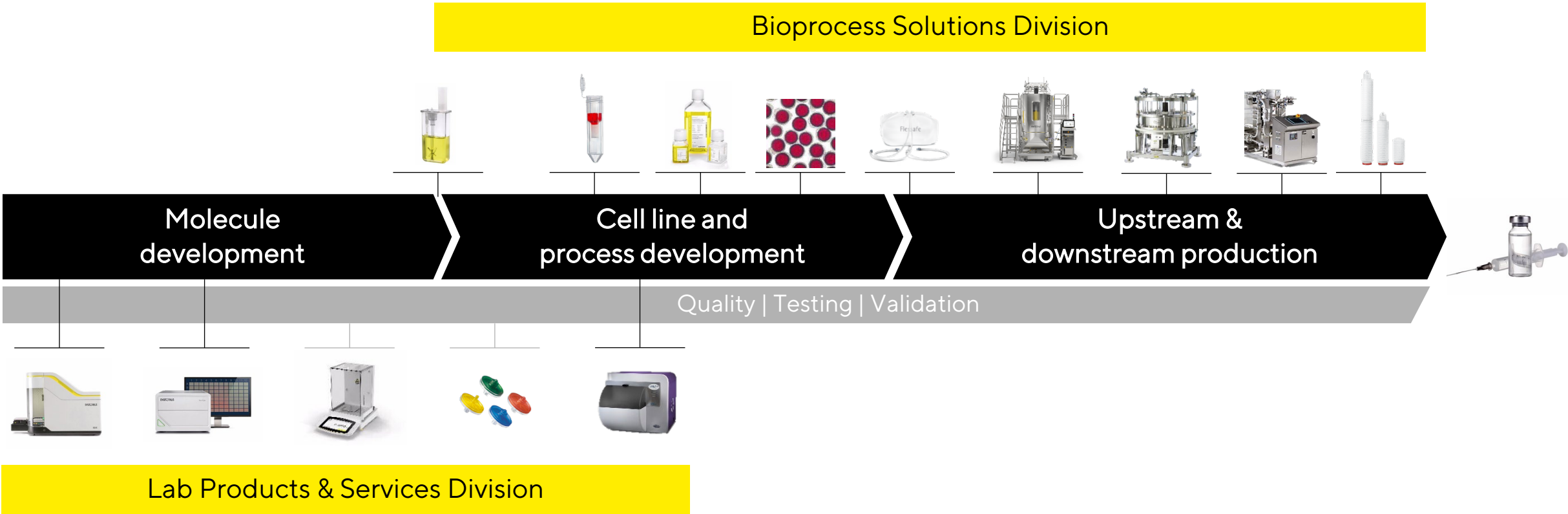
 **29.6%**
EBITDA margin^{1,2}

 **~€23.6bn**
Sartorius AG market capitalization;
included in relevant indices in Germany



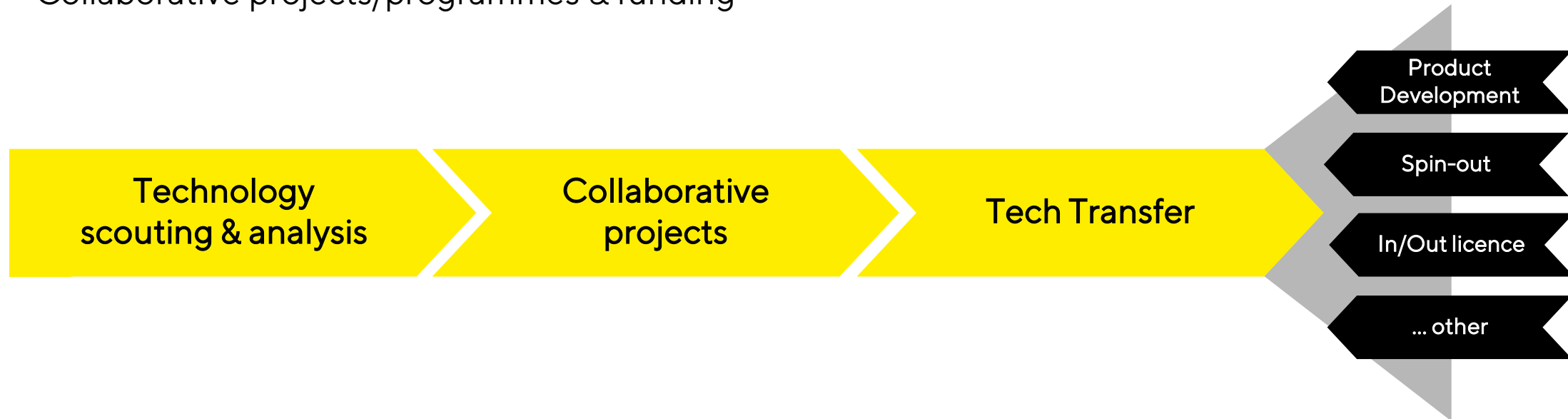
1 As of December 31, 2020, 2 Underlying EBITDA

Strategic focus on the biopharma market



Sartorius Corporate Research

- New ideas to concepts & novel applications aligned with Sartorius strategy (mission/vision)
- Securing of early stage IP positions
- Identify, nurture & enable technology creation with an understanding of end use
- Network with leading academia, consortia, research industry partner and opinion leaders
- Collaborative projects/programmes & funding



Summary

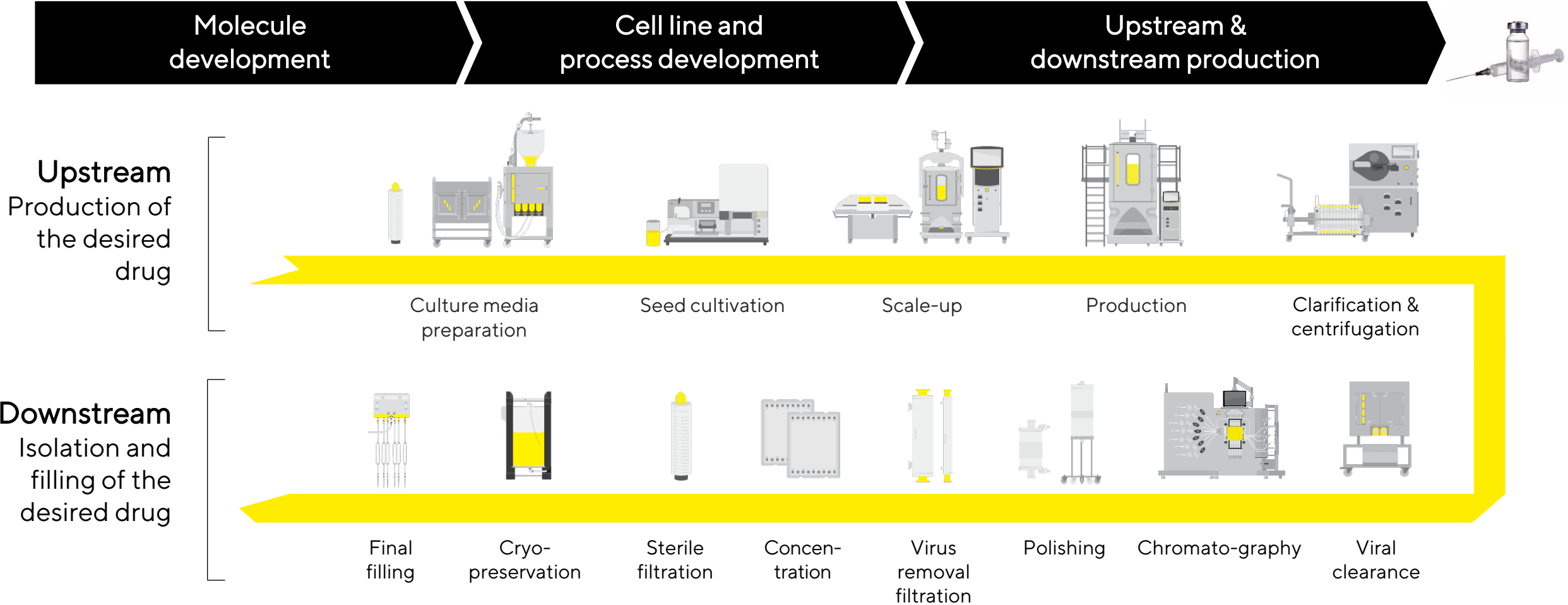
WHAT	WHY	HOW
Current practices in upstream bioprocessing	Shifting paradigms in biopharma R&D strategy	Making value out of data using modeling tools
The Biosimulator a generic upstream hybrid modeling structure	Unifying workflows from product development to manufacturing	Future perspectives and reflections on in-silico bioprocessing

WHAT	WHY	HOW
Current practices in upstream bioprocessing	Shifting paradigms in biopharma R&D strategy	Making value out of data using modeling tools
The Biosimulator a generic upstream hybrid modeling structure	Unifying workflows from product development to manufacturing	Future perspectives and reflections on in-silico bioprocessing

Manufacturing of biopharmaceuticals

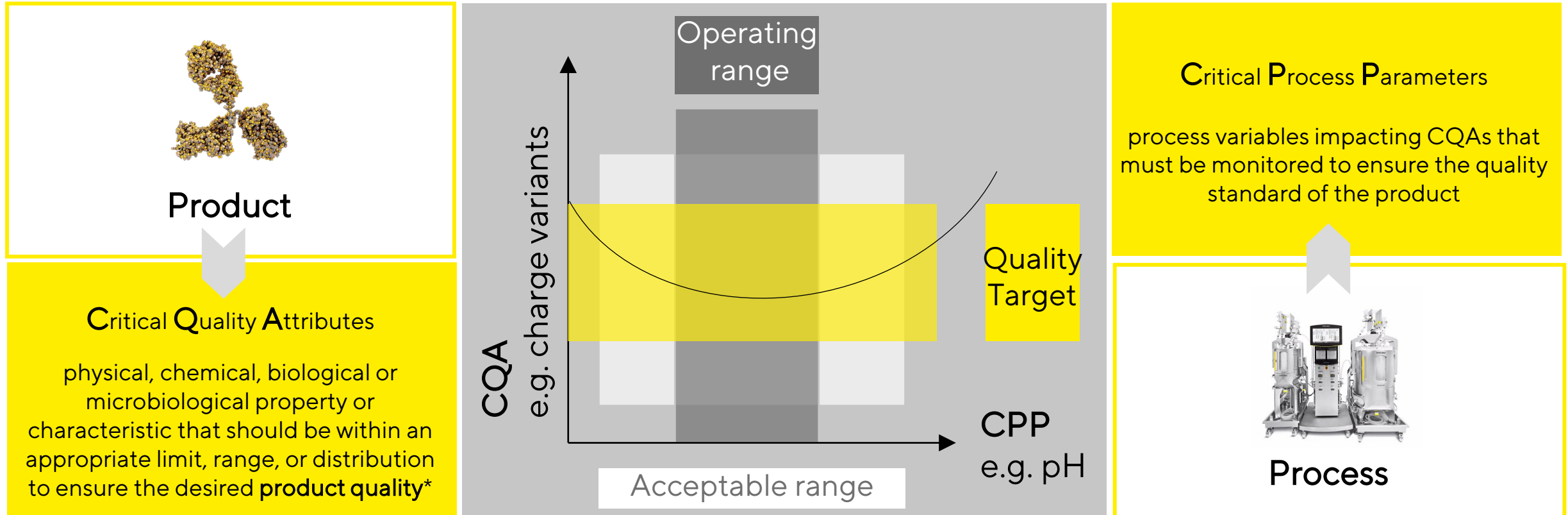


Large molecules
(> 20k atoms)
produced by
living cells culture



Bioprocess Performance Evaluation

CPPs are determined to ensure consistent and reproducible CQAs acquisition



While multiple researches have highlighted the potential benefit to shift the focus **from CPPs to CQAs**, process performance is still mainly evaluated in **process space** using Risk Analysis, DOE and MVDA

WHAT	WHY	HOW
Current practices in upstream bioprocessing	Shifting paradigms in biopharma R&D strategy	Making value out of data using modeling tools
The Biosimulator a generic upstream hybrid modeling structure	Unifying workflows from product development to manufacturing	Future perspectives and reflections on in-silico bioprocessing

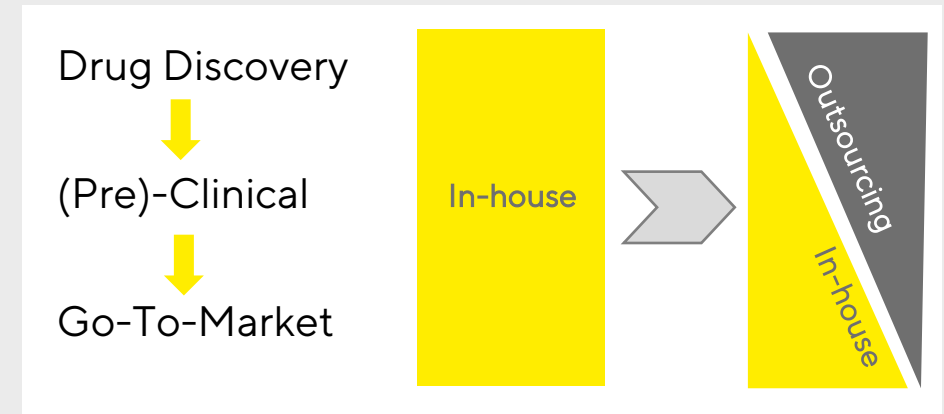
Challenging the biopharma R&D strategy

R&D cost are growing faster over an average investment period than actual revenues over the same period

This **NOT** due to a lack of Innovation or Market

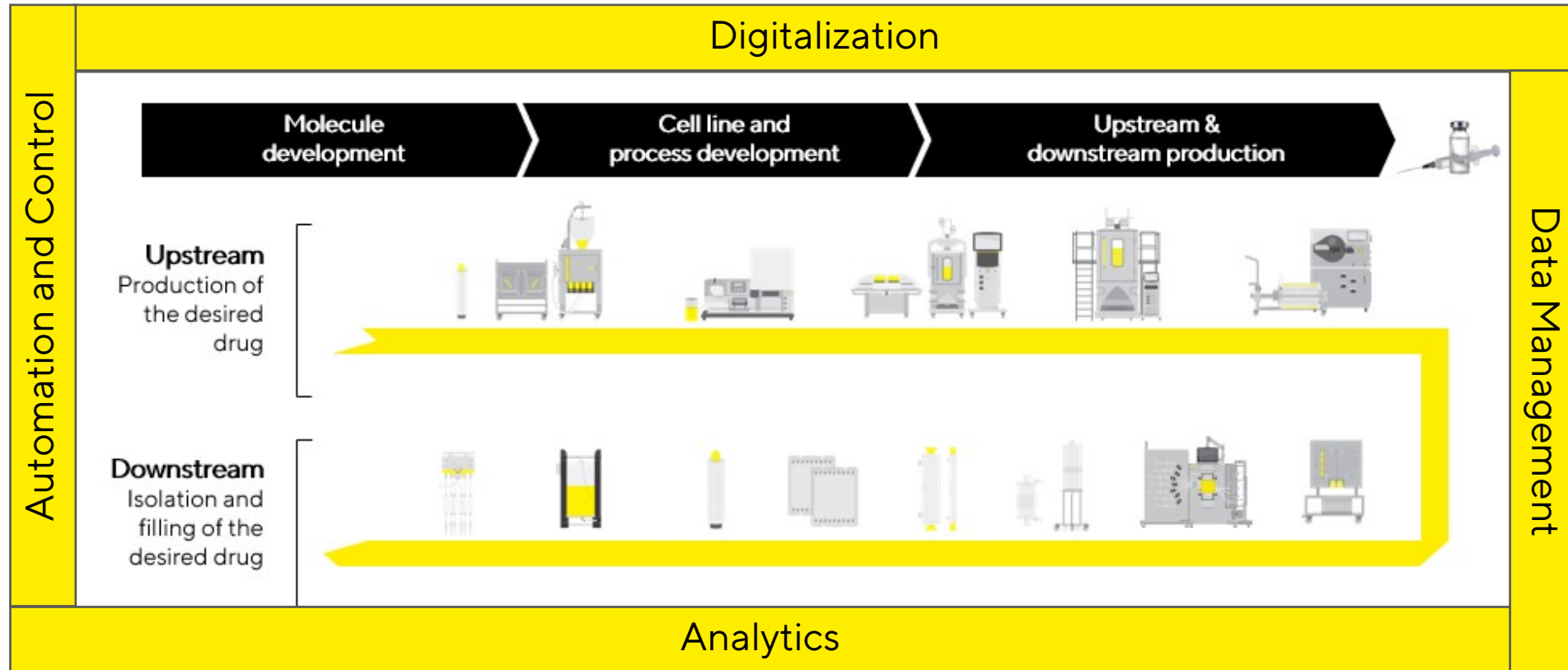
Law of diminishing returns ? Kelvin Stott

Inherent unpredictability of drug discovery and outdated business model? Standish Fleming



Outsourcing to external entrepreneurial and research community to lower the cost and increase the agility to face market disruption

Acceleration of Operational Tasks



WHAT	WHY	HOW
Current practices in upstream bioprocessing	Shifting paradigms in biopharma R&D strategy	Making value out of data using modeling tools
The Biosimulator a generic upstream hybrid modeling structure	Unifying workflows from product development to manufacturing	Future perspectives and reflections on in-silico bioprocessing

Process DOT* in the biopharmaceutical industry

Status

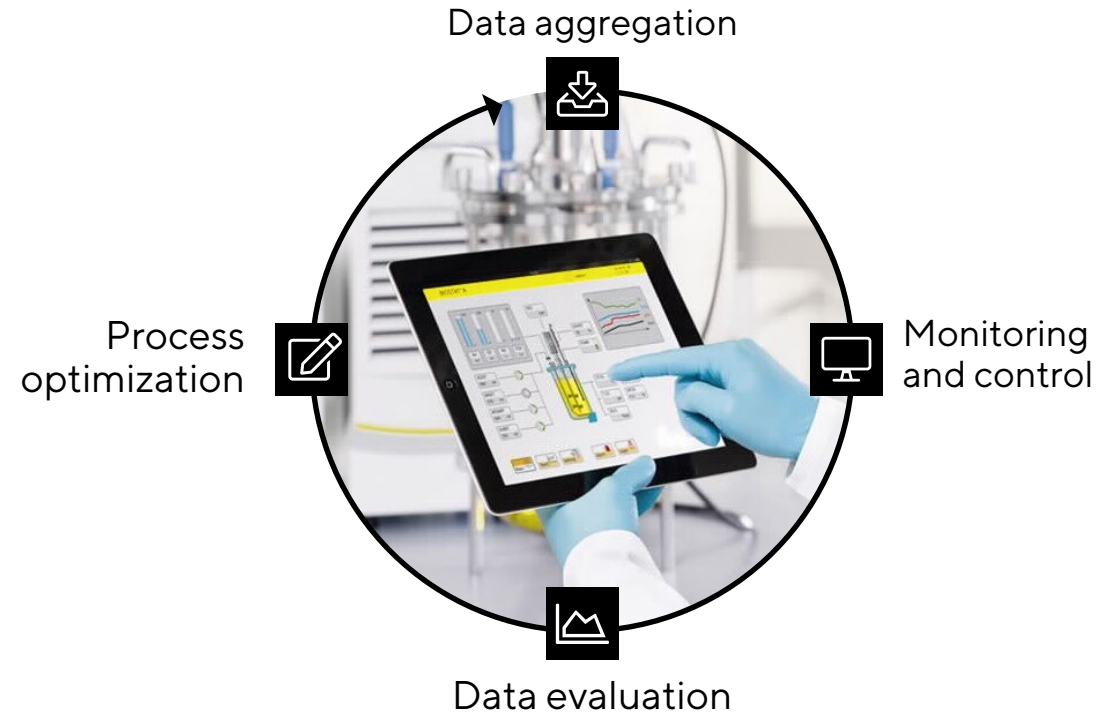
- Largely trial and error development
- Large amounts of complex heterogeneous data
- Long, costly development times

Challenges

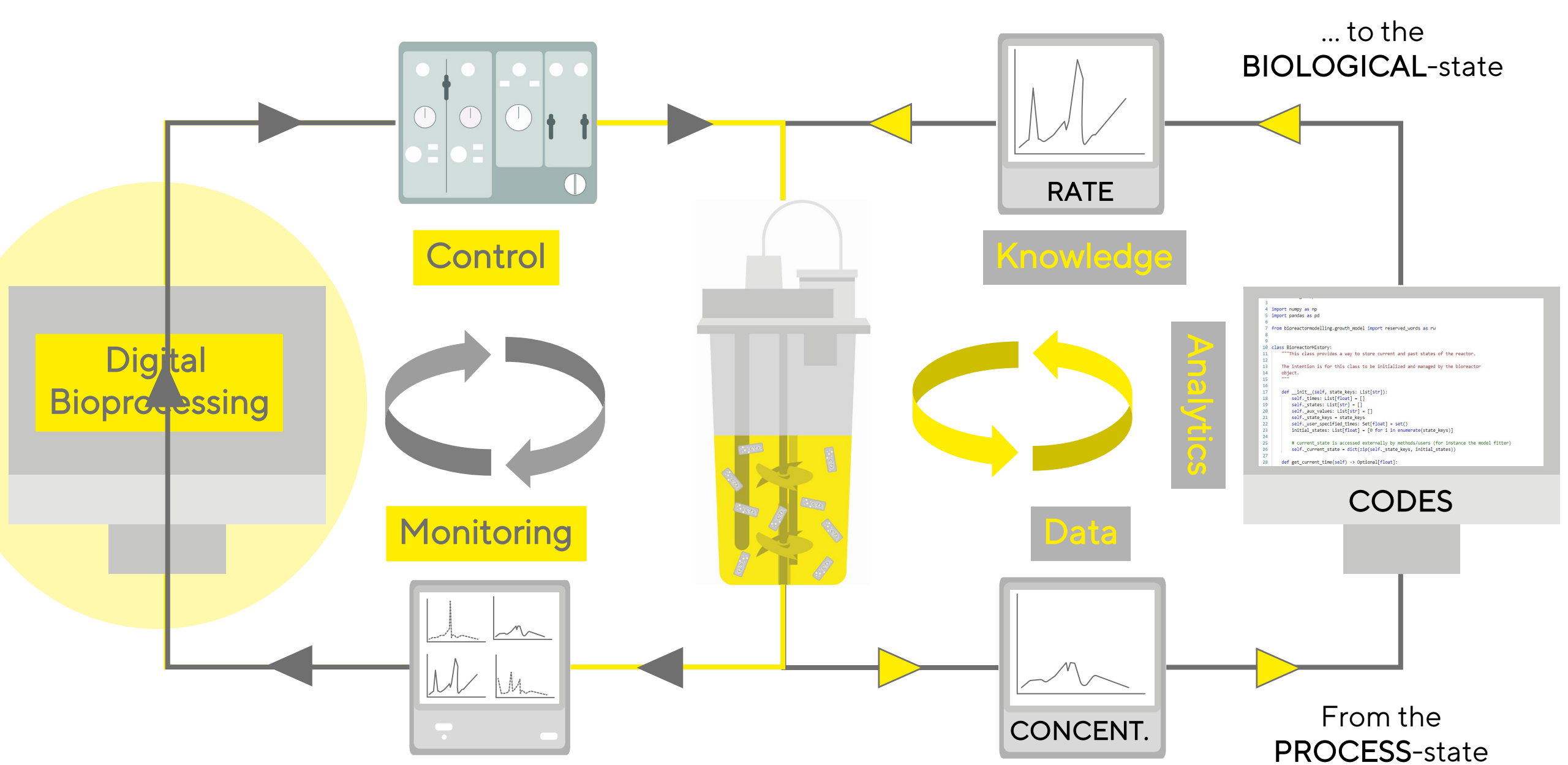
- Complex biology not fully understood
- Non-stationary, non-linear and stochastic behaviors
- Standard workflows do not yet exist

Opportunities

- In-silico modeling and simulation
- Real-time analytics and process optimization



- ✓ Rapid prototyping and process development
- ✓ Improved process performance and product quality
- ✓ De-risked transfer to manufacturing



WHAT	WHY	HOW
Current practices in upstream bioprocessing	Shifting paradigms in biopharma R&D strategy	Making value out of data using modeling tools
The Biosimulator a generic upstream hybrid modeling structure	Unifying workflows from product development to manufacturing	Future perspectives and reflections on in-silico bioprocessing

The Biosimulator

Substrate



Model calibrated from measured data
Transferrable to new cell lines

Hybrid State Space

Mechanistic
Process

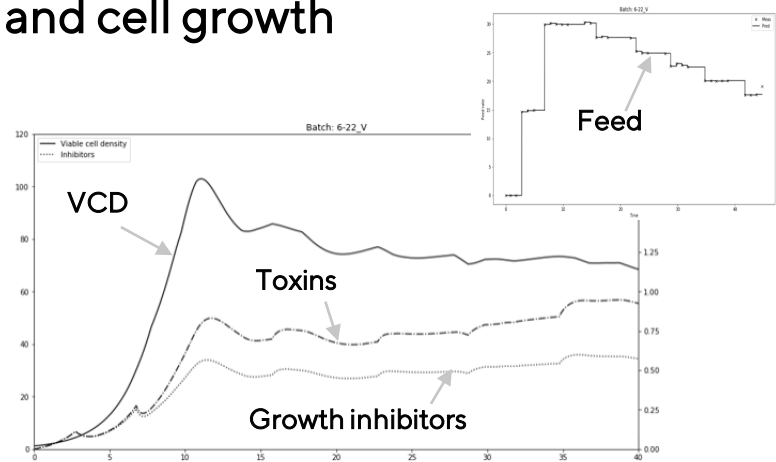
Data-driven
Metabolic



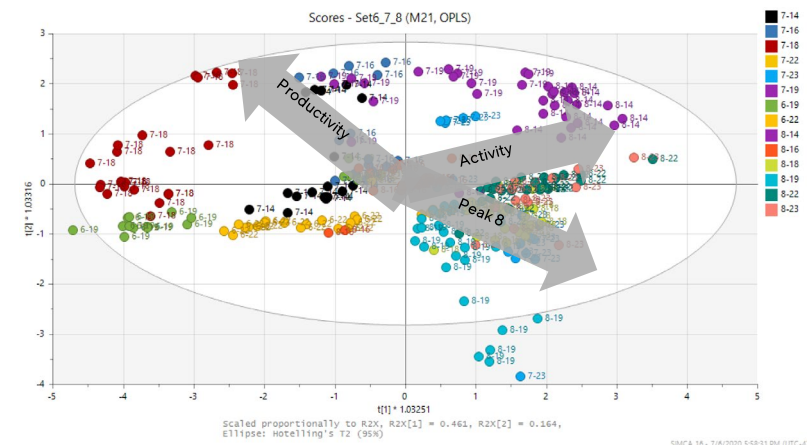
Product



Link between process operation
and cell growth



What are the cells doing?

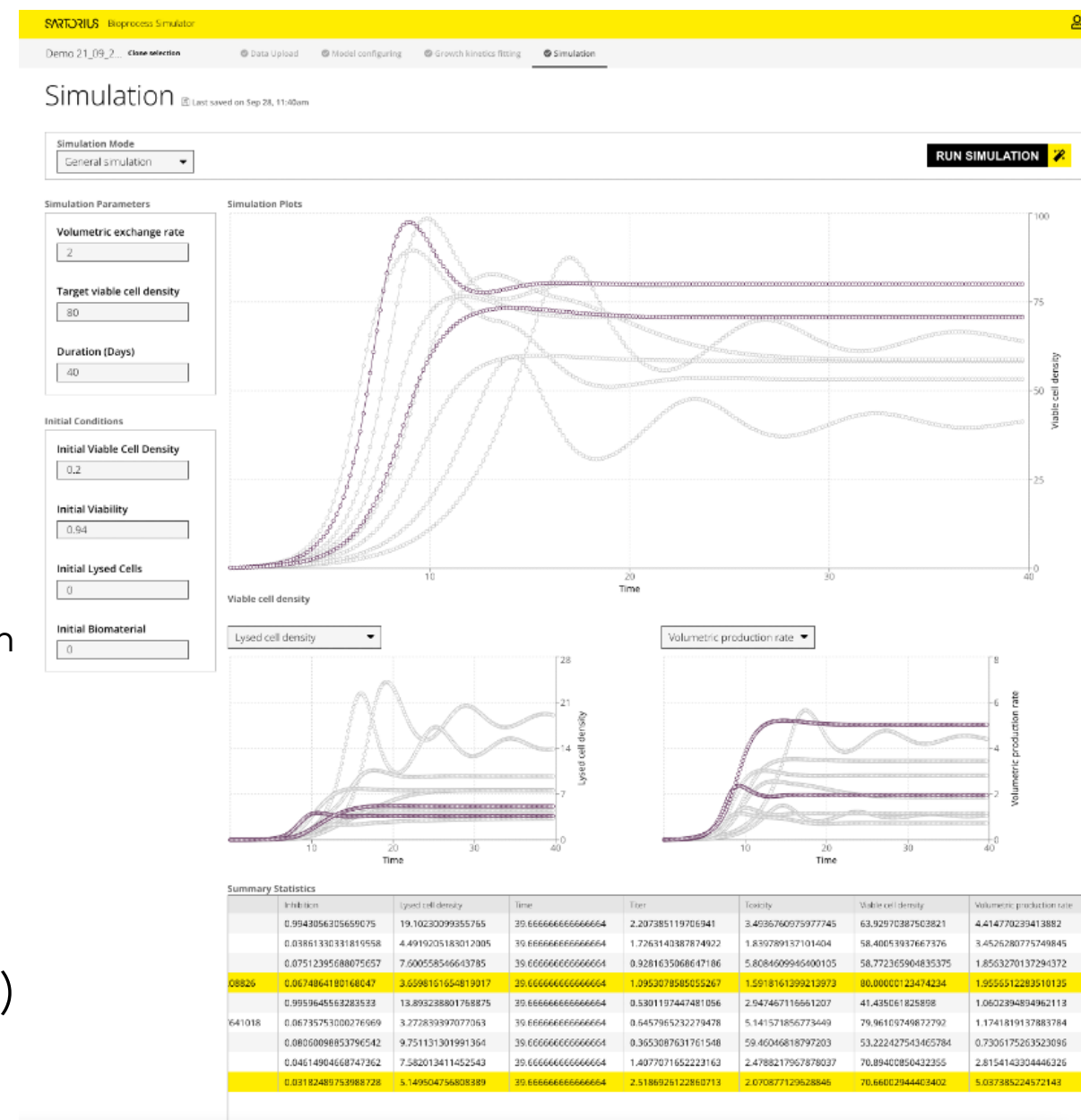


Commercial Product for Process Development

■ Two workflows

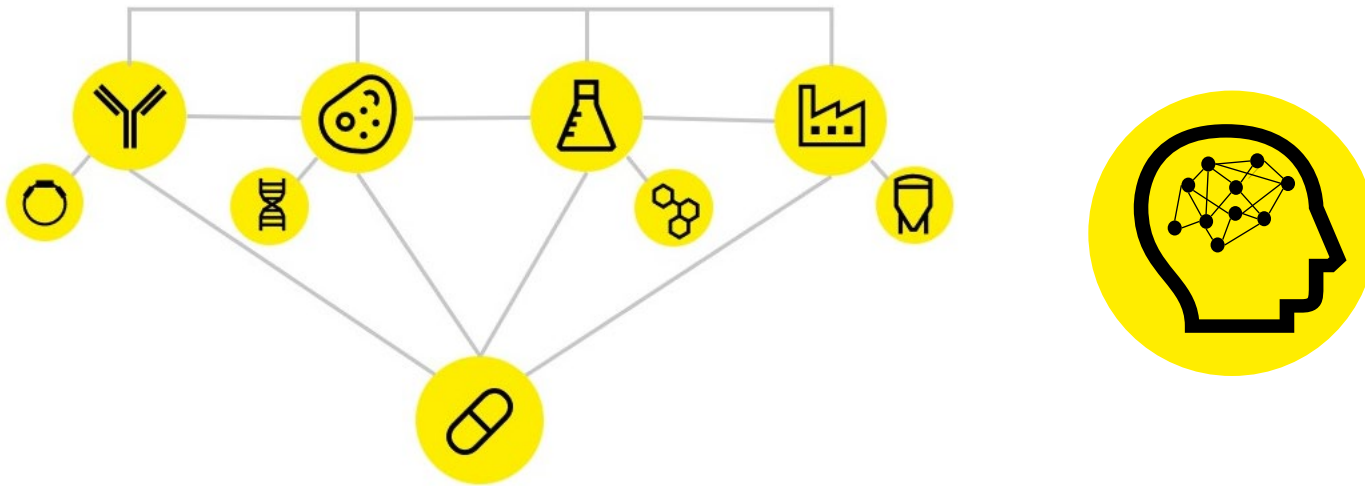
- ❑ Clone selection for intensified and perfusion
 - Forecast multiple cell line's performance in media exchange
- ❑ Digital twin for *in-silico* experimentation
 - What-if analysis: test impact of process adjustments on growth and product yield
 - Operating mode: transfer between fed-batch, intensified and perfusion
 - Media optimization
 - Critical Quality Attribute predictions

- Web/cloud based application (PC/tablet/iPad)
 - Integration in Umetrics Studio (2022)



WHAT	WHY	HOW
Current practices in upstream bioprocessing	Shifting paradigms in biopharma R&D strategy	Making value out of data using modeling tools
The Biosimulator a generic upstream hybrid modeling structure	Unifying workflows from product development to manufacturing	Future perspectives and reflections on in-silico bioprocessing

In-silico Experimentation Workflow



Process Development

- Enhanced clone selection
- Digital Twin: *Insilico* experimentation
- Optimization (fed-batch or intensified/continuous)

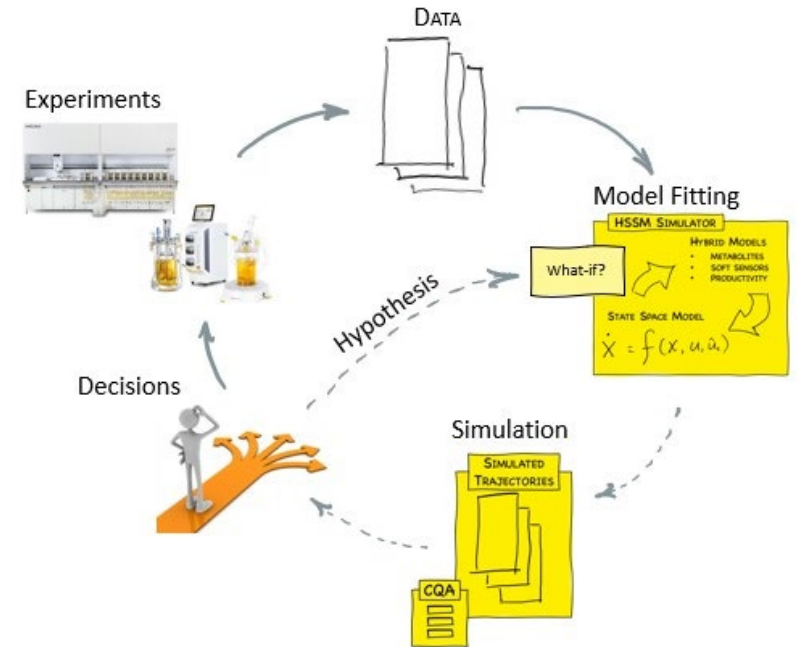
Manufacturing / Advanced control

- Intensified/continuous control
- Soft sensors
- Supervisory control (up/down-stream)

Engineering Studies

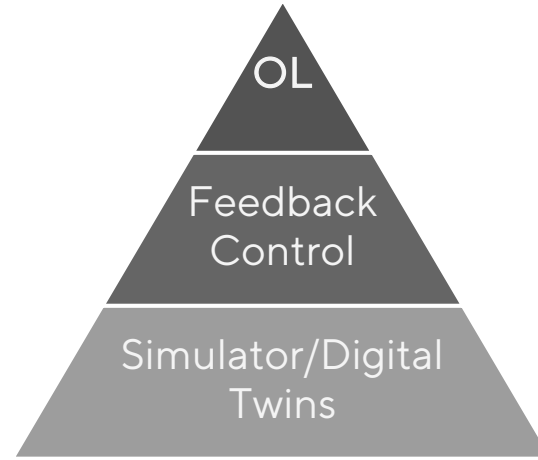
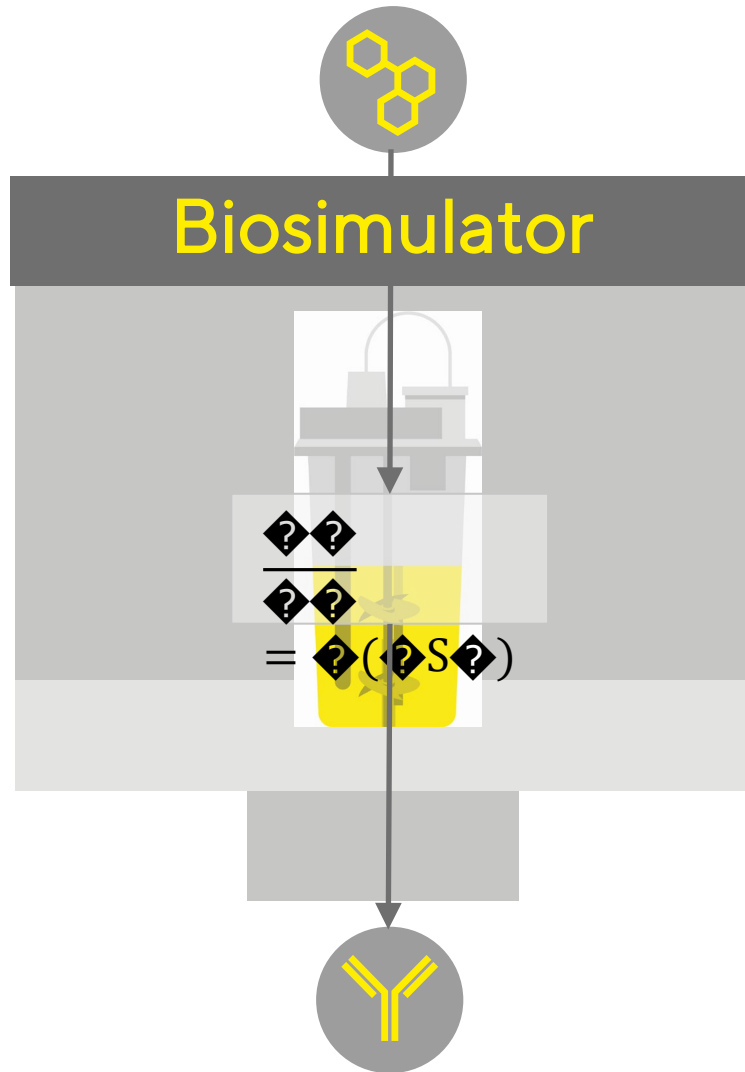
- Sizing and configuration of unit operations
- Plant costing
- End-to-end economic optimization

Systems thinking

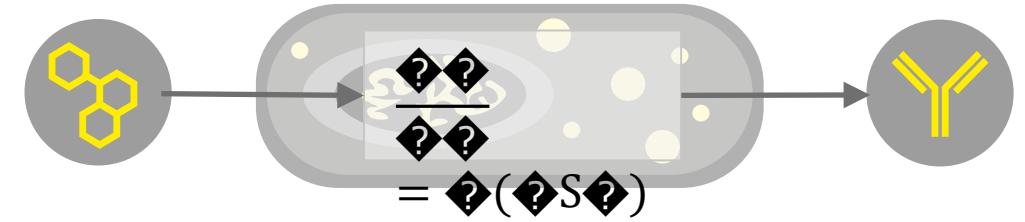


WHAT	WHY	HOW
Current practices in upstream bioprocessing	Shifting paradigms in biopharma R&D strategy	Making value out of data using modeling tools
The Biosimulator a generic upstream hybrid modeling structure	Unifying workflows from product development to manufacturing	Future perspectives and reflections on in-silico bioprocessing

Research Roadmap



Biological Systems modeling



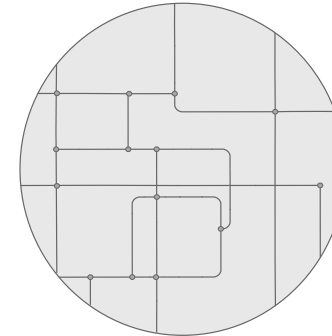
Process development

Analyze of process parameters influence



Macro

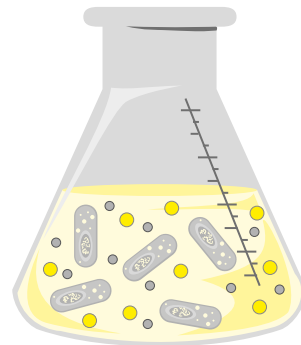
Optimization and control



Process engineering

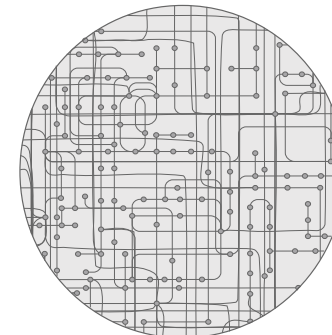
Media design

Identification of potential growth booster



Micro

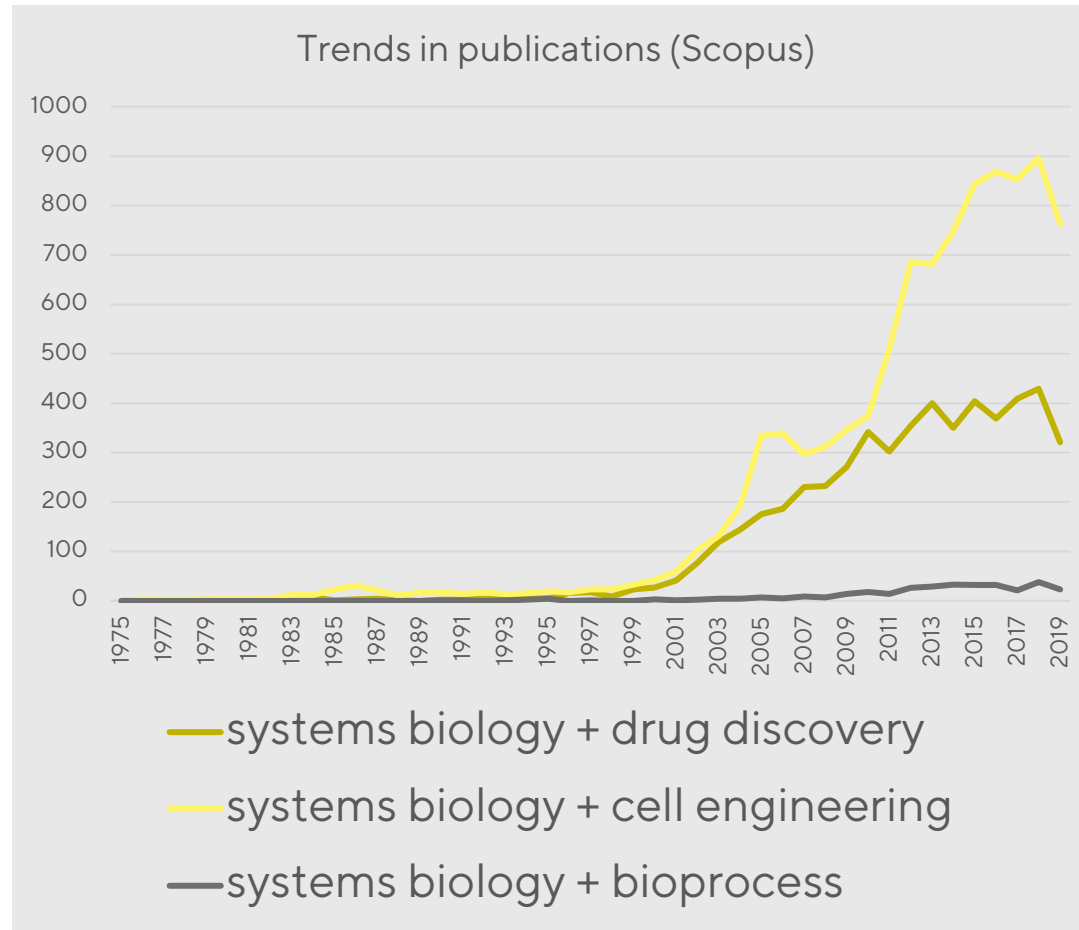
Description and understanding



Systems Biology

Systems Biology for Process DOT*

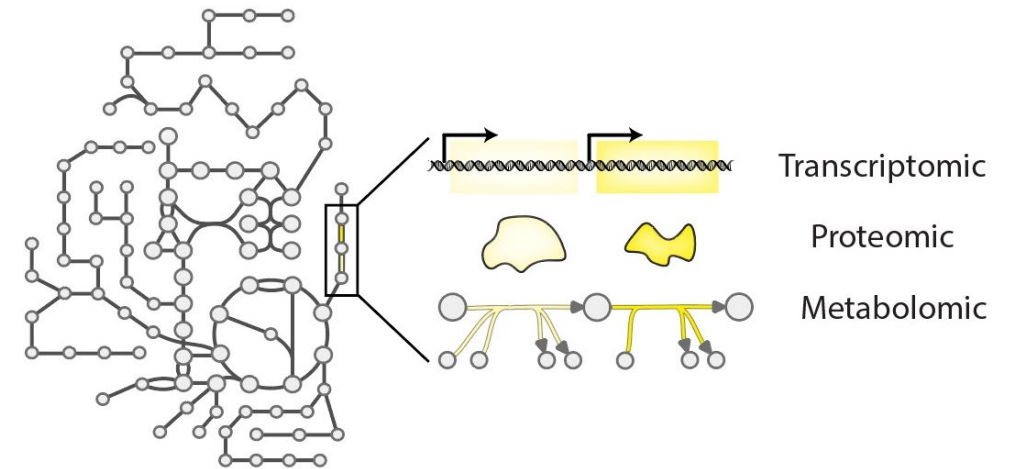
Note : Multiple reasons have been highlighted as main factors currently limiting the adoption of systems biology tools**



Metabolic Networks

enable the coherent organization of large datasets into biological networks

→ provide non-intuitive insights on biological systems that *in vivo* experiments alone cannot provide



→ link between genotype and phenotype

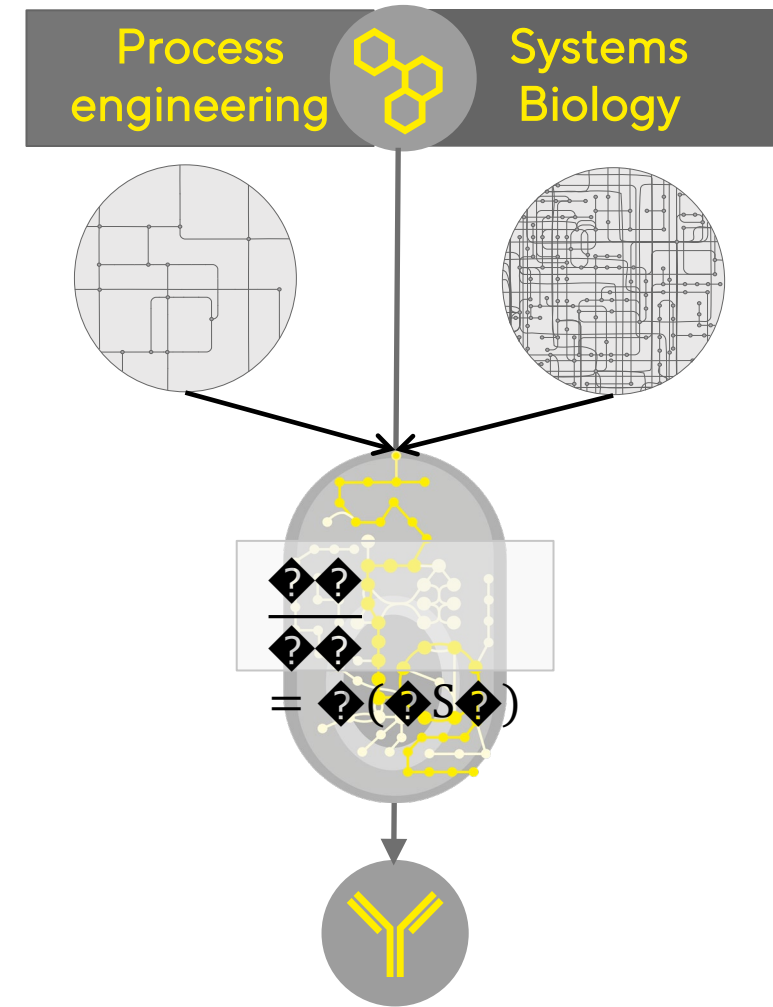
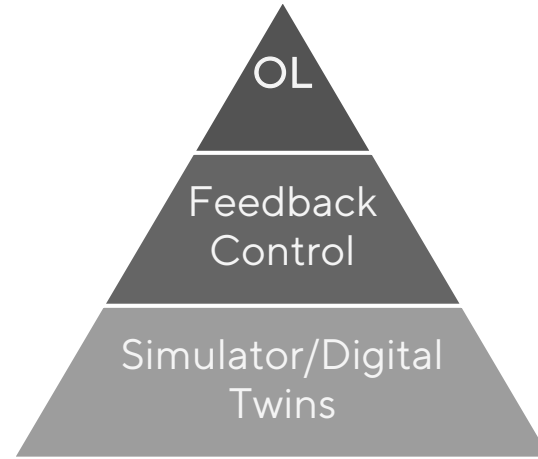
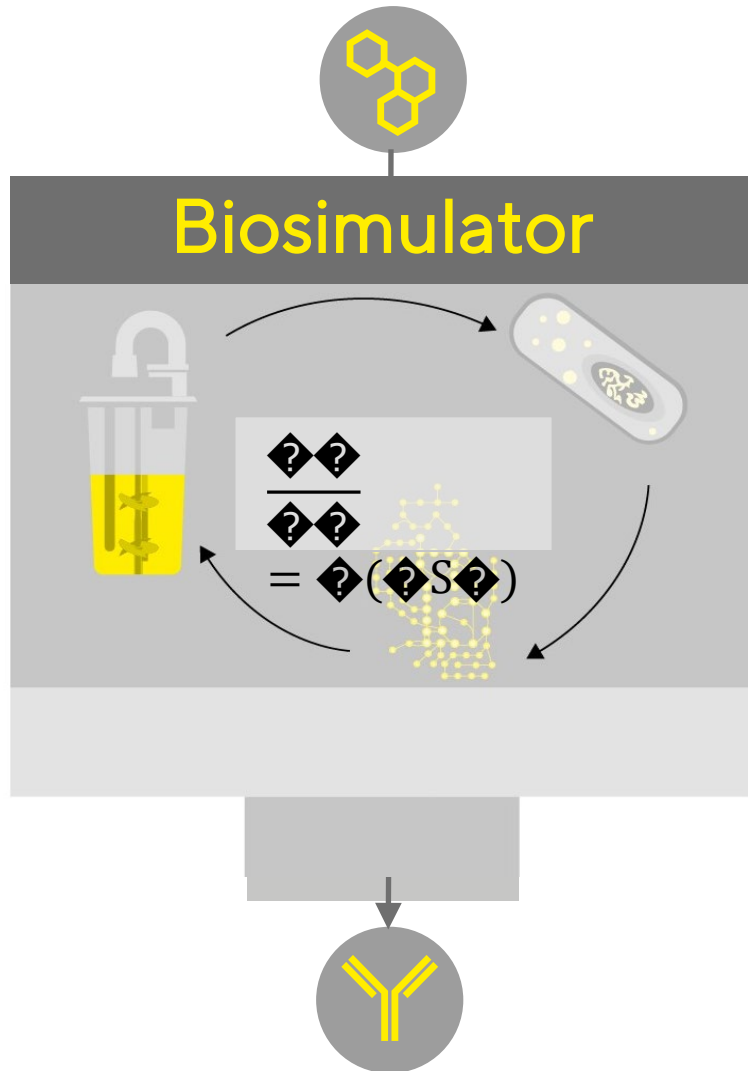
*DOT = Design, Optimization and Transfer

** Richelle et al. *npj Syst Biol Appl* 6, 6 (2020).

Towards a widespread adoption of metabolic modeling tools in biopharmaceutical industry: a process systems biology engineering perspective.

SARTORIUS

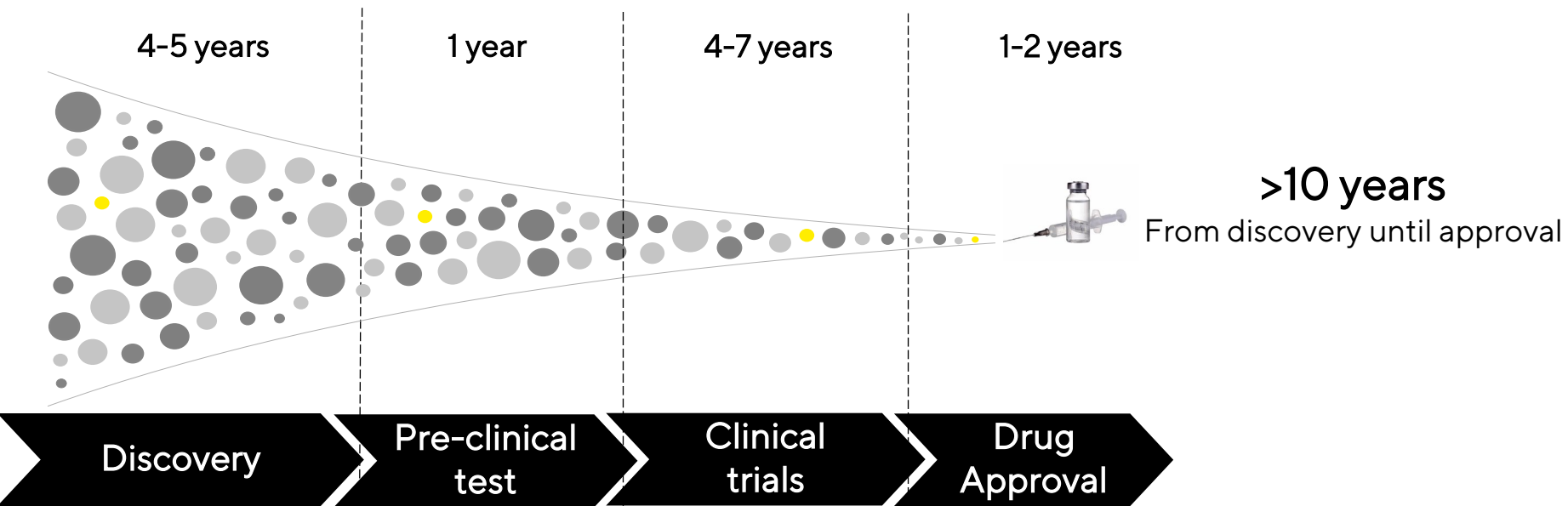
Research Roadmap



WHAT	WHY	HOW
Current practices in upstream bioprocessing	Shifting paradigms in biopharma R&D strategy	Making value out of data using modeling tools
The Biosimulator a generic upstream hybrid modeling structure	Unifying workflows from product development to manufacturing	Future perspectives and reflections on in-silico bioprocessing

Future outlook on drug development

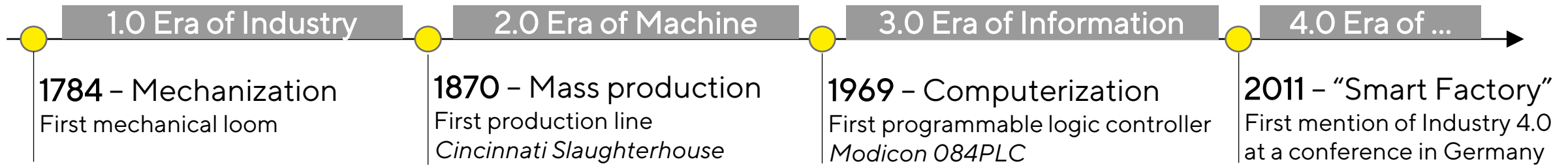
Only one out of 10,000 new drug candidates reaches the market



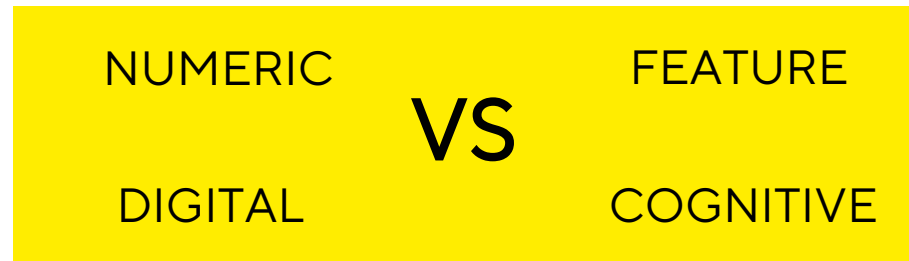
>€2bn
Average costs of developing
a successful drug

~10%
Probability of clinical success
(Phase I to approval)

Food for thoughts – Bioprocessing 4.0



“The [AI] field's heavy use of statistical techniques to pick regularities in masses of data is unlikely to yield the explanatory insight that science ought to offer”. Noam Chomsky, 2012



➔ Shifting existing paradigm for defining regulated biomanufacturing processes from data- to knowledge-based approaches

Thank you.

Acknowledgments

Advanced Data Analytics team

Olivier Cloarec
Brandon Corbett
Christoffer Edlund
Chris McCready
Emil Nylind
Rickard Sjoegren
Elsa Soerman Paulsson
Alexander Sutherland
Johan Trygg
Anton Verneresson

Anne Richelle
Senior Research Scientist
Sartorius Corporate Research
anne.richelle@sartorius.com

SARTORIUS