

# SARTORIUS

## Simplifying Progress

### Putting the “D” in CDMO with Data Analytics and Ambr<sup>®</sup>

Tiffany McLeod

Kevin McHugh



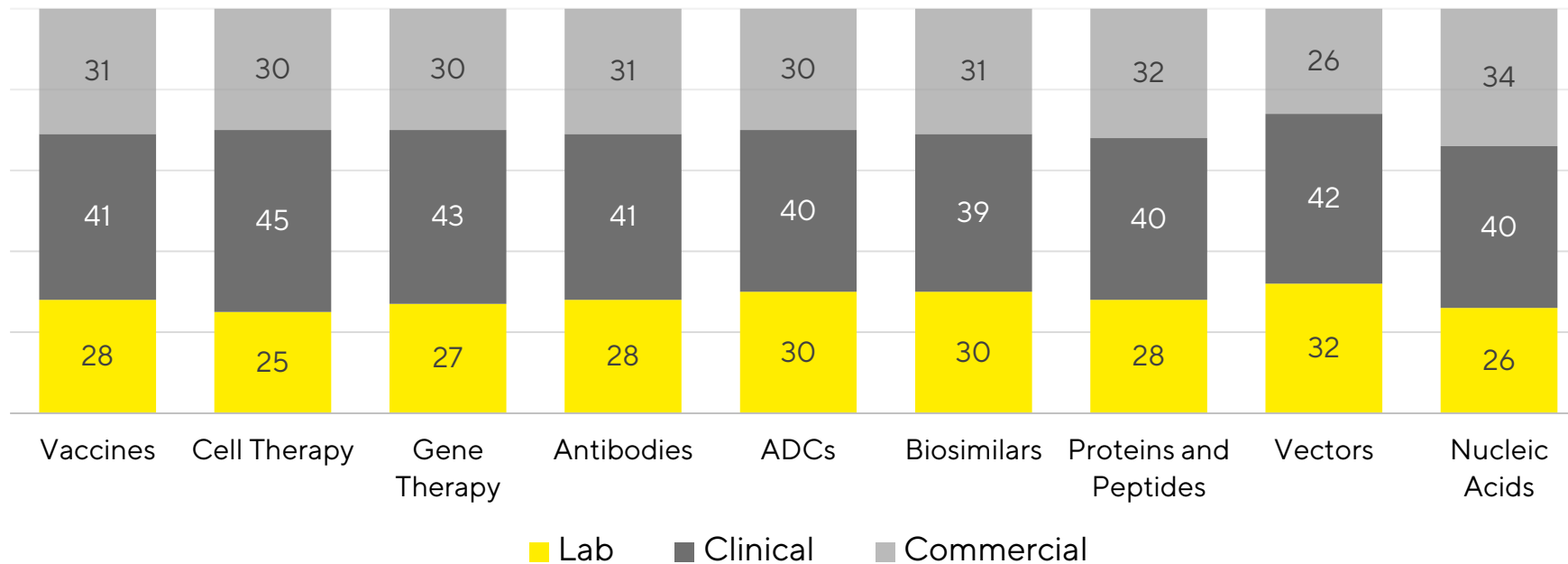
# Agenda

- Market Trends & Analysis
- CDMOs Pursuing QbD
- Technologies that Can Accelerate Process Development at CDMOs
  - Ambr<sup>®</sup> Systems Overview
- Why Ambr<sup>®</sup> and Data Analytics?



# Biopharmaceutical CDMOs: Distribution of Biologics by Operations

## INTER-SCALE DISTRIBUTIONS



- Lab and Clinical operations make up a majority of the CDMO operations
- 57% of lab scale, 56% of clinical scale operations have been occupied by vaccines, antibodies, and proteins and peptides

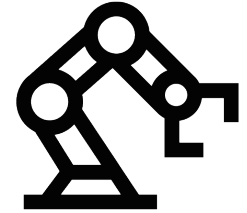
Roots Analysis, Biopharmaceutical Contract Manufacturing Market (3<sup>rd</sup> Edition) 2019-2030 (2019)

# Bioprocess Development Technology Trends

**In-Silico Experimentation** This disruptive technology influencing all industries including Pharma and Biotech and the hype is real. In-silico-based tools are part of drug targeting, screening and discovery, clinical studies and predictive analytics related to any risk involved.



**High-throughput Process Development** Involves the miniaturization, automation, and parallelization of process development activities – proving for a systematic approach for time- and resource-efficient workflow.



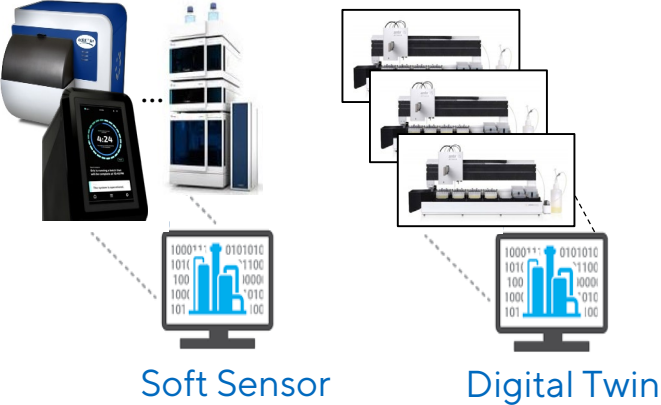
**Continuous Bioprocessing** The intensification of both upstream and downstream operations will require higher levels of control during PD and presents new scale-up challenges.





# Bioprocess Development Technology Trends

## In-Silico Experimentation



## High-throughput Process Development



## Continuous Bioprocessing



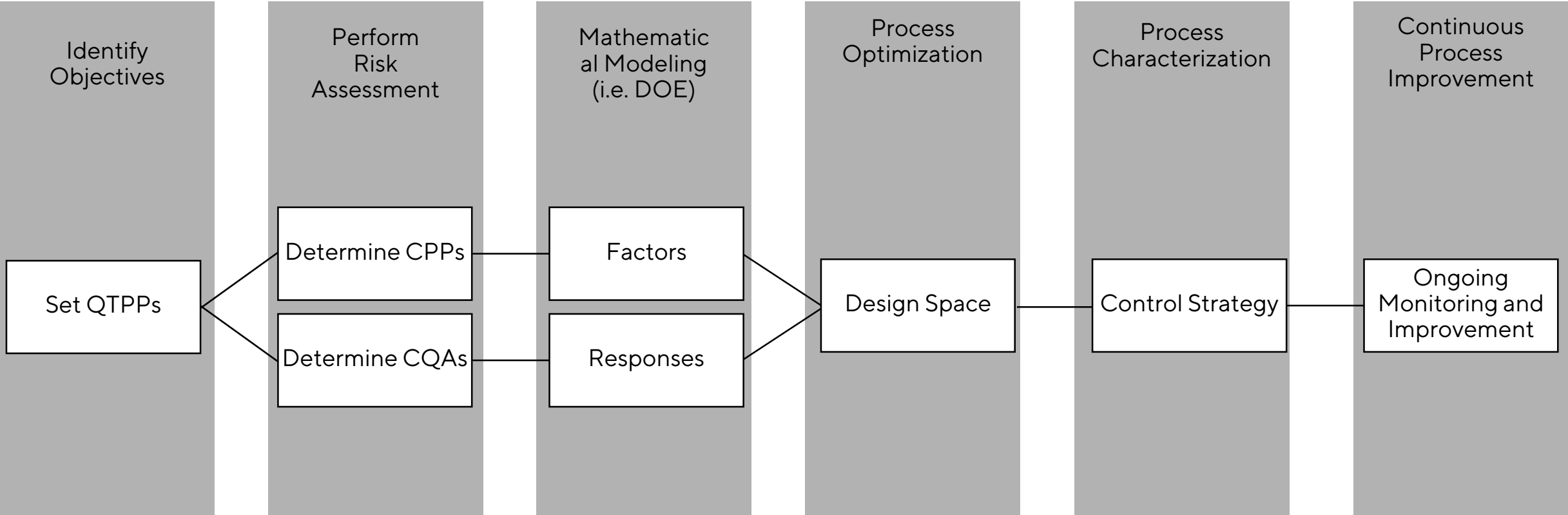
# Agenda

- Market Trends & Analysis
- CDMOs Pursuing QbD
- Technologies that Can Accelerate Process Development at CDMOs
  - Ambr<sup>®</sup> Systems Overview
- Why Ambr<sup>®</sup> and Data Analytics?





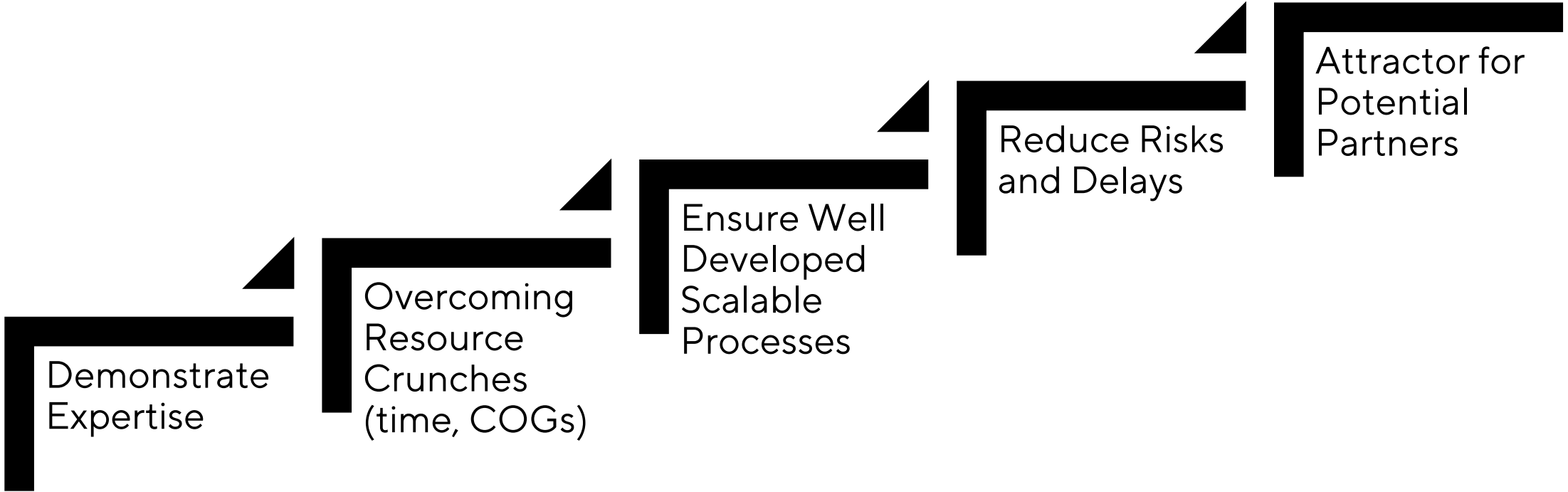
# Why Regulatory Agencies are Pushing Quality by Design (QbD)



"QbD means designing and developing manufacturing processes during the product development stage to consistently ensure a predefined quality at the end of the manufacturing process." ICH Q10, FDA 2006



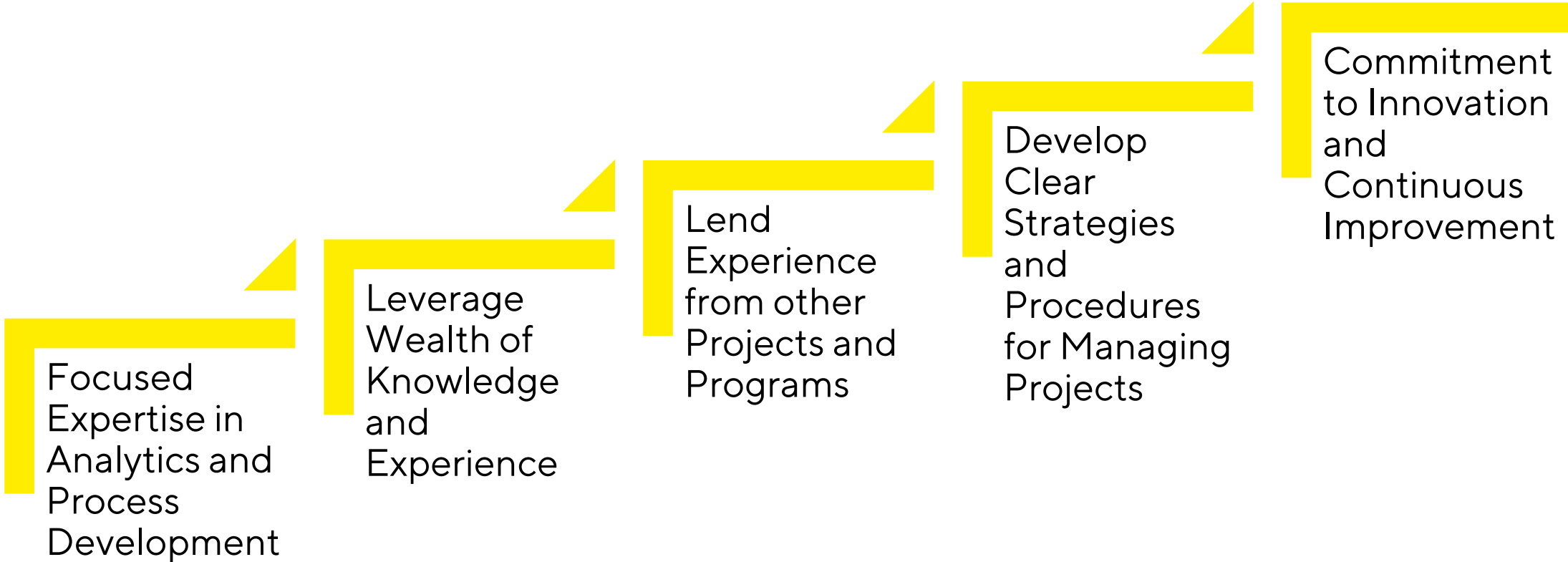
# Why Should CDMOs Have Strong QbD Packages?







# How Can CDMOs Develop Strong QbD Packages?



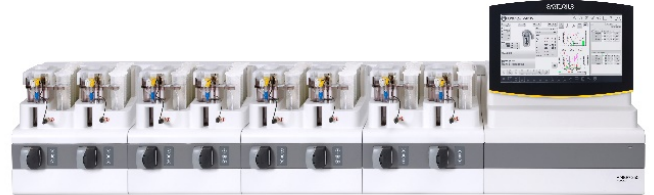
# Agenda

- Market Trends & Analysis
- CDMOs Pursuing QbD
- Technologies that Can Accelerate Process Development at CDMOs
  - Ambr<sup>®</sup> Systems Overview
- Why Ambr<sup>®</sup> and Data Analytics?





# Fully Scalable Range of Single-Use Bioreactors from 15 mL to 2000 L Following Conventional Stirred Tank Design Principles





# Ambr<sup>®</sup> 15 Cell Culture



More representative →



← Predictive



- Clone selection
- Media and feed screening
- Early process optimization
- Batch, fed-batch and perfusion mimic



- Transient transfection studies
- Mammalian and insect cell lines
- Suspension cell lines and adherent cell lines on microcarriers (e.g 293, Vero)
- Applications ranging from recombinant protein vaccines to viral vectors and mRNA

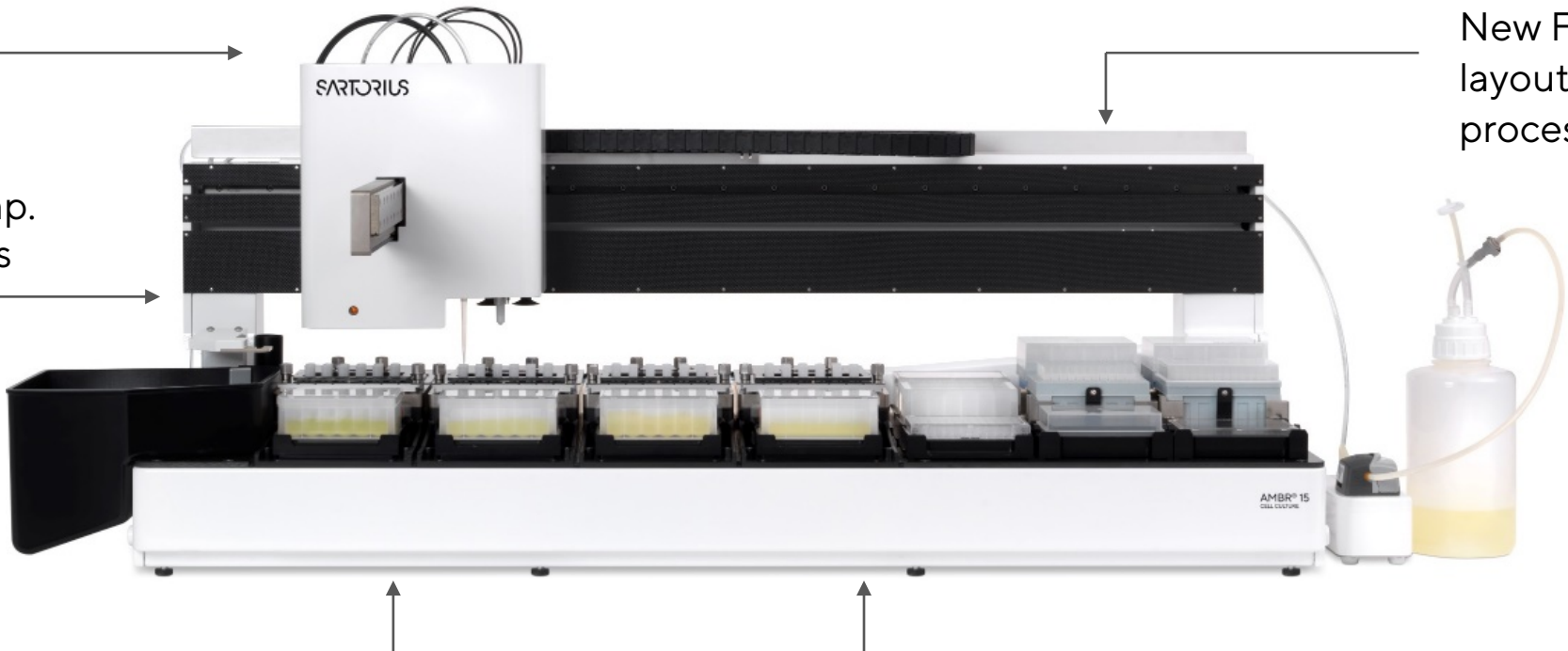


# Ambr<sup>®</sup> 15 Generation 2 Workstation Features and Functionality

New Liquid Handler improves precision

New Culture Station design improves temp. control and increases stirrer speed range

New Flexible deck layout allows more process flexibility



New Large tip bin expands tip capacity

24 or 48 vessels arranged in 2 or 4 culture stations

Active cooling option

New Rapid vessel drain allows removal of culture or spent media from microbioreactors



# Enabling Simple Consistent Clone/Strain Selection



Currently the selection of the top clones or strains is performed upon visual inspection of specific quality parameters or one number readout (e.g. max titre or IVCC)

The selection process is user dependent and the full power of using all available data with a balanced tool to give a robust priority selection is **not fully utilized**

Achieve a **better** and **more robust** clone selection in a **standardized** way using a balanced priority of CQA's



# Ambr<sup>®</sup> 250 High Throughput: the Fast Track to Intensified Cell Culture Process Development

This next-generation bioreactor system can enable and enhance your perfusion process development in these key areas:

- Reduce complexity and setup time with single-use perfusion bioreactors
- Increase experiment capacity and data consistency with a fully automated parallel system
- Gain predictive, scalable results with the industry standard Ambr<sup>®</sup> 250 bioreactor
- Increase cost effectiveness due to media and labor cost savings
- Improve development timelines and performance with larger studies and perfusion DoE



# Ambr<sup>®</sup> 250 High Throughput Workstations

12 Vessel

24 Vessel

Perfusion Option

Analysis Module

Automated system for aseptic bolus addition and sampling

13 feet

Class II equivalent BSC



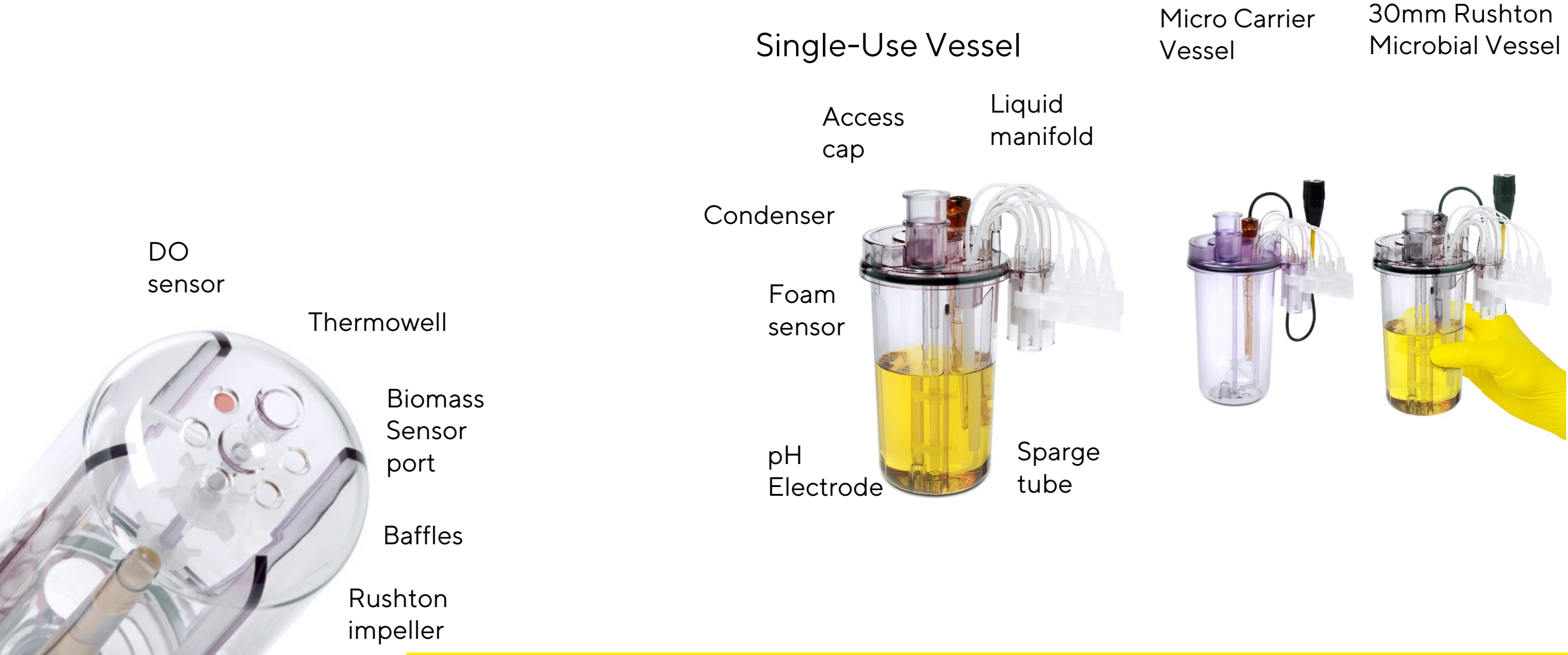
Chilled sample storage with an option for freezer positions

Bed locations for pipette tips and stirred bed location for 1 L and 175 mL bottles

12 or 24 bioreactors 4 pumps each

Perfusion pump towers and filters

# Ambr<sup>®</sup> 250 High Throughput Vessels



# Ambr<sup>®</sup> 250 High Throughput Perfusion Option– Key Features

- Fits into the existing footprint of a standard ambr<sup>®</sup>250 high throughput system
- Fully capable of operating in either fed-batch or perfusion modes
- ATF & TFF consumable includes a single-use perfusion filter and single-use pump chambers
- Includes a high efficiency sparger for enhanced oxygen transfer and capable of supporting high cell densities
- Single-use bleed and permeate bags
- Medium exchange of 0.5 - 4 VVD
- Integrated automation





# Ambr<sup>®</sup> 250 Modular System

## Single-Use Vessel

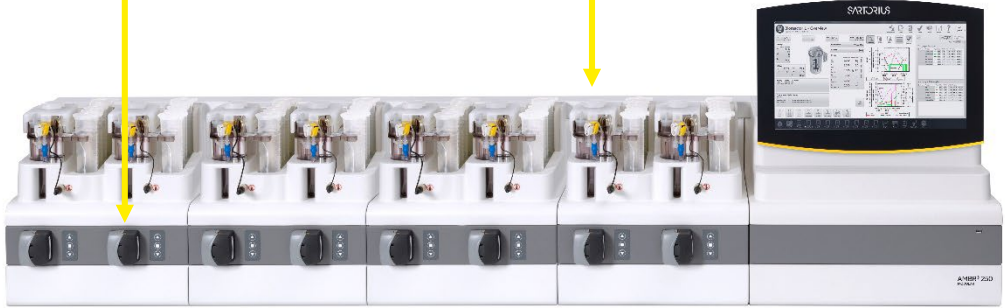


Integrated off-gas analyzers available for CO<sub>2</sub>, and O<sub>2</sub> Measurement, as well as OUR, CER, and RQ calculation

## Workstation

Variable speed 'x pump' allows for flexible use and quick reactor harvesting.

Simple and efficient pH and gassing, and agitation connections



# Easily Develop Advanced Control Strategies

- Flexible Control Strategies:
  - Setpoints can be based on a value, equation, profile, or even OPC input.
  - Not just for pumps but also process parameters
- Easy to use Triggers
  - Variety of options available to easily automate the transition between phases of a process
- Flexible Control Loops:
  - Any process parameter or calculated value can have a setpoint and an output can be assigned to control it
  - Implement advanced control strategies like DO-Stat in just a few clicks
- Calculated Variables
  - Equations can be developed to calculated values based on various sensors and inputs
  - Can be used as triggers, controlled variables & PID terms

The screenshot displays a complex software interface for developing advanced control strategies. It features several overlapping windows:

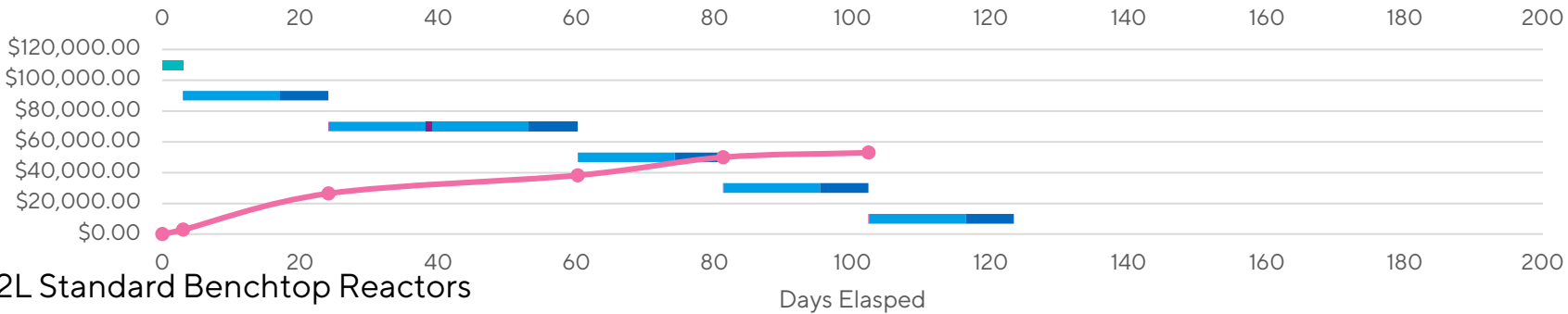
- Edit Expression:** Shows an expression  $OUR / (CalcKla\_Cstar - CalcKla\_C\_Broth)$  and a status "Expression is valid". It includes a variable selection grid with categories like "Standard operators" and "Variables".
- Edit step - 'Start Pump Feed#1':** Contains the text "Edit step parameters" and "Start pump running and set the profile for the pump".
- Edit custom variable 'CalcKla\_kla':** Shows "Edit parameters for variable" and a table for defining a PID loop.
- Edit loop 'kLa' controlling CalcKla\_kla.SP:** Displays "Parameters for a set of cascaded PID loops" with a table of properties and values.

Property	Value
Description	
Description	kLa
Level 1 - Stir speed	
Output	Stir speed
Effect of output	Increases controlled variable
Dead band	
Minimum (rpm)	800
Maximum (rpm)	4500
Proportional Term - kP	10
Derivative Time - tD (s)	0
Integral Time - tI (s)	100

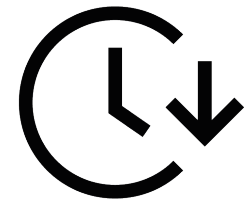
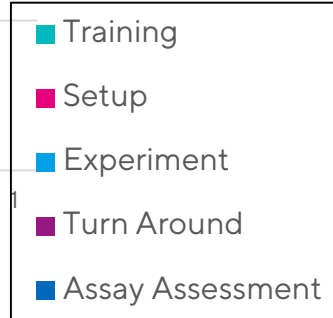
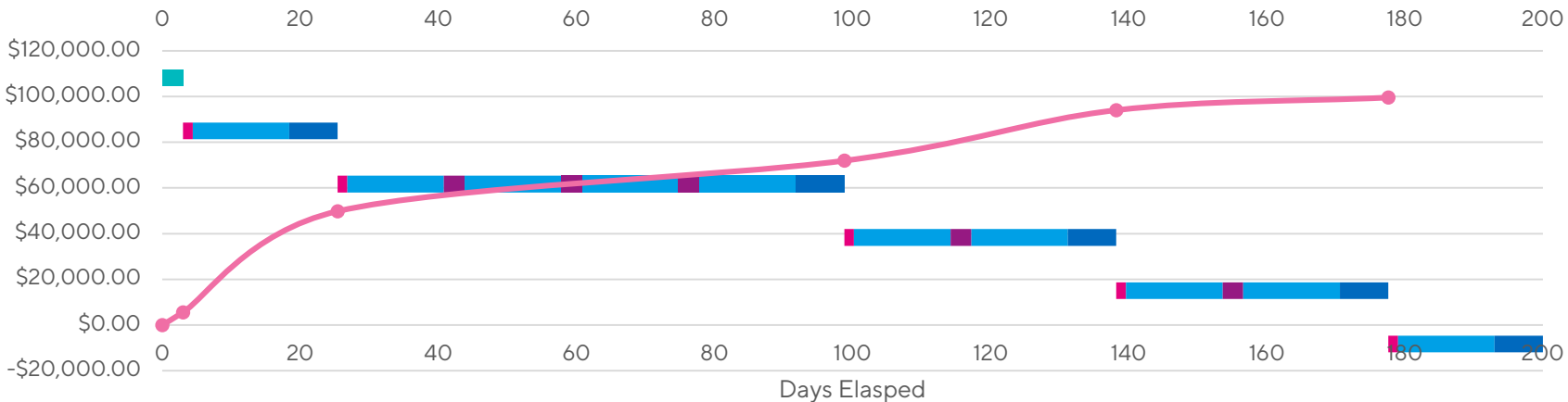


# High Throughput Upstream Process Characterization Platform Development – Cost and Time Saving Comparison With Traditional Approach

## ■ Ambr® 250 – 24 Vessel Configuration



## ■ 12x2L Standard Benchtop Reactors



# Scalable, Automated At-Line and Real Time Process Monitoring and Control

From 15 mL to 2000 L SU bioreactors

Guidance for Industry  
PAT — A Framework for  
Innovative Pharmaceutical  
Development, Manufacturing,  
and Quality Assurance

PAT facilitates the  
implementation of  
QbD!



NOVA Flex II  
with ESM



ViCell XR

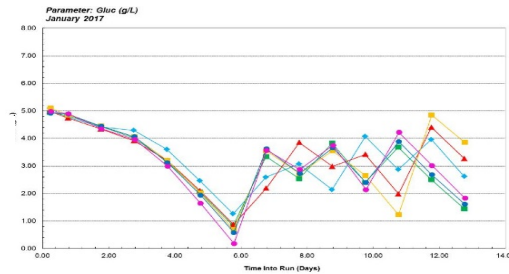
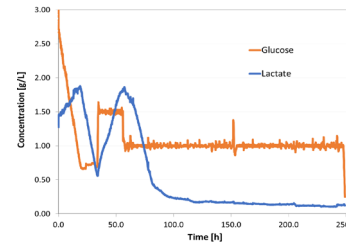


Off-gas  
and Foam  
Sensors

Online  
Biomass



Analysis  
Module  
w/ BioPAT  
Spectro



BioPAT® Trace



BioPAT® X-gas



BioPAT® Foam



EC/SU pH

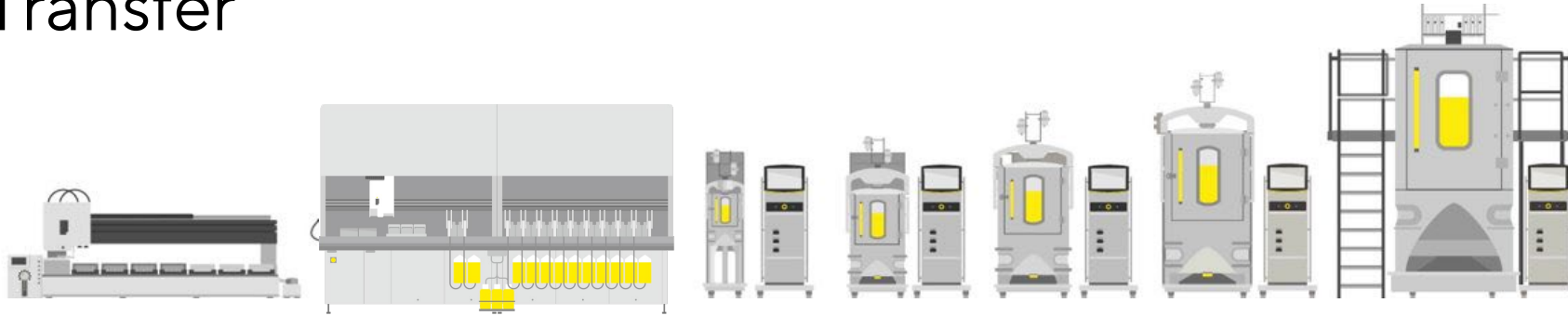


BioPAT® Viamass

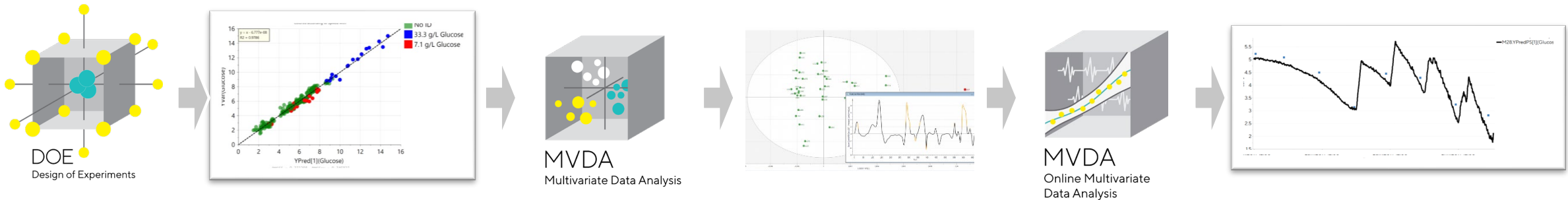


BioPAT® Spectro

# Spectroscopy Platform across Scales enabling straight-forward Model-Transfer



Rowland-Jones, RC, Graf, A, Woodhams, A, et al. Spectroscopy integration to miniature bioreactors and large scale production bioreactors—Increasing current capabilities and model transfer. *Biotechnol Progress*. 2020;e3074. <https://doi.org/10.1002/btpr.3074>



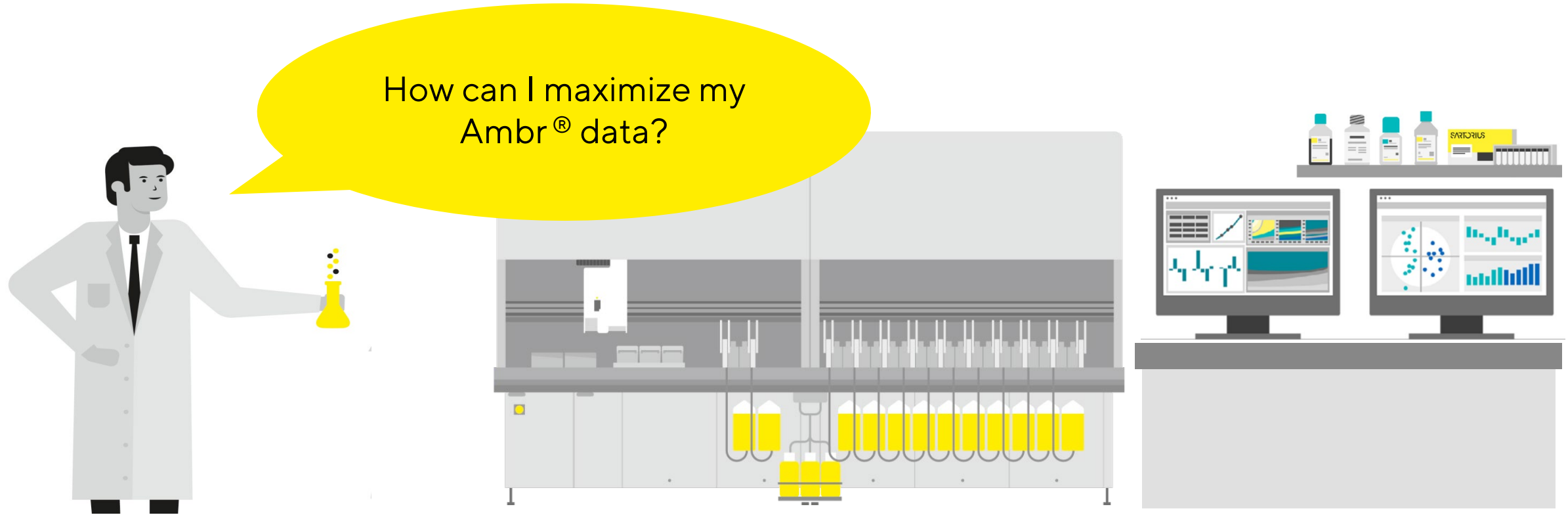


# Agenda

- Market Trends & Analysis
- CDMOs Pursuing QbD
- Technologies that Can Accelerate Process Development at CDMOs
  - Ambr<sup>®</sup> Systems Overview
- Why Ambr<sup>®</sup> and Data Analytics?



# Ambr<sup>®</sup> and Data Analytics: The Perfect Match



# From Data to Decisions

Sartorius Data Analytics tools examine large amounts of data to uncover hidden patterns.

MODDE®



SIMCA®



# Design, Execute and Analyze Experiments Quickly and Easily

- MODDE® Design of Experiments
  - Familiar interface for defining factors and responses.
  - Functionality to automatically import setpoints and extract responses from ambr15 experiments
    - Not only end point but min, max, timepoints, etc.
  - Software automatically accounts for system type (12, 24, 48 & culture station limitations)
- Experiments and Results Viewer
  - Quickly and easily explore previous experiments and campaigns
  - Quickly generate tables and charts
  - Investigate experiment audit trails
  - Combine and visualize data from multiple experiments
- Easy to Use SIMCA® Data Export
  - Quickly and easily export online and at-line sequential batch data for an experiment or entire campaign
  - Interpolate data as needed

The screenshot displays the SIMCA software interface. The 'Factor Definition' window is open, showing 'Factor name' and 'Units' fields. The 'Results/Bioreactors 1-2' window shows a line graph of 'red pixels' vs 'time' for 'Bioreactor selection'. An 'Export plotted data' dialog box is overlaid, allowing users to export data for the lines plotted on the graph. The dialog includes options for 'Data points' (All), 'Export format' (Stack bioreactors (for SIMCA)), 'Phase column' (None), 'Date format' (Local times), 'Interpolation' (Interpolated), 'Interpolation method' (Mean value), and 'Interpolation interval' (5m). Buttons for 'Save settings...', 'Load settings...', 'Ok', and 'Cancel' are visible. In the background, a 'Results/Tables' window shows a table of data for various bioreactors.

Bioreactor	Time (hour)	Value
Bioreactor 1	1.51	1.51
Bioreactor 1	30.01	30.01
Bioreactor 1	43.78	43.78
Bioreactor 1	53.00	53.00
Bioreactor 1	67.80	67.80
Bioreactor 1	77.96	77.96
Bioreactor 1	89.98	89.98
Bioreactor 1	91.80	91.80
Bioreactor 1	100.01	100.01
Bioreactor 1	107.99	107.99
Bioreactor 1	115.78	115.78
Bioreactor 1	124.03	124.03
Bioreactor 1	131.99	131.99
Bioreactor 1	139.85	139.85
Bioreactor 1	147.82	147.82
Bioreactor 1	155.99	155.99
Bioreactor 1	163.71	163.71
Bioreactor 1	172.03	172.03
Bioreactor 1	180.50	180.50
Bioreactor 1	190.01	23
Bioreactor 1	203.78	2.7
Bioreactor 1	211.97	3.1
Bioreactor 1	216.82	3.9
Bioreactor 1	227.78	5.0
Bioreactor 1	236.03	5.4
Bioreactor 1	244.01	6.4
Bioreactor 1	251.89	7.3
Bioreactor 1	259.87	7.7
Bioreactor 1	269.56	7.97
Bioreactor 1	278.01	8.1
Bioreactor 1	286.22	8.2
Bioreactor 1	294.29	8.3
Bioreactor 1	302.19	8.4
Bioreactor 1	310.00	8.5
Bioreactor 1	317.68	8.6
Bioreactor 1	325.25	8.7
Bioreactor 1	332.71	8.8
Bioreactor 1	340.06	8.9
Bioreactor 1	347.30	9.0
Bioreactor 1	354.44	9.1
Bioreactor 1	361.48	9.2
Bioreactor 1	368.42	9.3
Bioreactor 1	375.26	9.4
Bioreactor 1	382.00	9.5
Bioreactor 1	388.64	9.6
Bioreactor 1	395.18	9.7
Bioreactor 1	401.62	9.8
Bioreactor 1	407.96	9.9
Bioreactor 1	414.20	10.0
Bioreactor 1	420.44	10.1
Bioreactor 1	426.68	10.2
Bioreactor 1	432.92	10.3
Bioreactor 1	439.16	10.4
Bioreactor 1	445.40	10.5
Bioreactor 1	451.64	10.6
Bioreactor 1	457.88	10.7
Bioreactor 1	464.12	10.8
Bioreactor 1	470.36	10.9
Bioreactor 1	476.60	11.0
Bioreactor 1	482.84	11.1
Bioreactor 1	489.08	11.2
Bioreactor 1	495.32	11.3
Bioreactor 1	501.56	11.4
Bioreactor 1	507.80	11.5
Bioreactor 1	514.04	11.6
Bioreactor 1	520.28	11.7
Bioreactor 1	526.52	11.8
Bioreactor 1	532.76	11.9
Bioreactor 1	539.00	12.0
Bioreactor 1	545.24	12.1
Bioreactor 1	551.48	12.2
Bioreactor 1	557.72	12.3
Bioreactor 1	563.96	12.4
Bioreactor 1	570.20	12.5
Bioreactor 1	576.44	12.6
Bioreactor 1	582.68	12.7
Bioreactor 1	588.92	12.8
Bioreactor 1	595.16	12.9
Bioreactor 1	601.40	13.0
Bioreactor 1	607.64	13.1
Bioreactor 1	613.88	13.2
Bioreactor 1	620.12	13.3
Bioreactor 1	626.36	13.4
Bioreactor 1	632.60	13.5
Bioreactor 1	638.84	13.6
Bioreactor 1	645.08	13.7
Bioreactor 1	651.32	13.8
Bioreactor 1	657.56	13.9
Bioreactor 1	663.80	14.0
Bioreactor 1	670.04	14.1
Bioreactor 1	676.28	14.2
Bioreactor 1	682.52	14.3
Bioreactor 1	688.76	14.4
Bioreactor 1	695.00	14.5
Bioreactor 1	701.24	14.6
Bioreactor 1	707.48	14.7
Bioreactor 1	713.72	14.8
Bioreactor 1	719.96	14.9
Bioreactor 1	726.20	15.0
Bioreactor 1	732.44	15.1
Bioreactor 1	738.68	15.2
Bioreactor 1	744.92	15.3
Bioreactor 1	751.16	15.4
Bioreactor 1	757.40	15.5
Bioreactor 1	763.64	15.6
Bioreactor 1	769.88	15.7
Bioreactor 1	776.12	15.8
Bioreactor 1	782.36	15.9
Bioreactor 1	788.60	16.0
Bioreactor 1	794.84	16.1
Bioreactor 1	801.08	16.2
Bioreactor 1	807.32	16.3
Bioreactor 1	813.56	16.4
Bioreactor 1	819.80	16.5
Bioreactor 1	826.04	16.6
Bioreactor 1	832.28	16.7
Bioreactor 1	838.52	16.8
Bioreactor 1	844.76	16.9
Bioreactor 1	851.00	17.0
Bioreactor 1	857.24	17.1
Bioreactor 1	863.48	17.2
Bioreactor 1	869.72	17.3
Bioreactor 1	875.96	17.4
Bioreactor 1	882.20	17.5
Bioreactor 1	888.44	17.6
Bioreactor 1	894.68	17.7
Bioreactor 1	900.92	17.8
Bioreactor 1	907.16	17.9
Bioreactor 1	913.40	18.0
Bioreactor 1	919.64	18.1
Bioreactor 1	925.88	18.2
Bioreactor 1	932.12	18.3
Bioreactor 1	938.36	18.4
Bioreactor 1	944.60	18.5
Bioreactor 1	950.84	18.6
Bioreactor 1	957.08	18.7
Bioreactor 1	963.32	18.8
Bioreactor 1	969.56	18.9
Bioreactor 1	975.80	19.0
Bioreactor 1	982.04	19.1
Bioreactor 1	988.28	19.2
Bioreactor 1	994.52	19.3
Bioreactor 1	1000.76	19.4
Bioreactor 1	1007.00	19.5
Bioreactor 1	1013.24	19.6
Bioreactor 1	1019.48	19.7
Bioreactor 1	1025.72	19.8
Bioreactor 1	1031.96	19.9
Bioreactor 1	1038.20	20.0
Bioreactor 1	1044.44	20.1
Bioreactor 1	1050.68	20.2
Bioreactor 1	1056.92	20.3
Bioreactor 1	1063.16	20.4
Bioreactor 1	1069.40	20.5
Bioreactor 1	1075.64	20.6
Bioreactor 1	1081.88	20.7
Bioreactor 1	1088.12	20.8
Bioreactor 1	1094.36	20.9
Bioreactor 1	1100.60	21.0
Bioreactor 1	1106.84	21.1
Bioreactor 1	1113.08	21.2
Bioreactor 1	1119.32	21.3
Bioreactor 1	1125.56	21.4
Bioreactor 1	1131.80	21.5
Bioreactor 1	1138.04	21.6
Bioreactor 1	1144.28	21.7
Bioreactor 1	1150.52	21.8
Bioreactor 1	1156.76	21.9
Bioreactor 1	1163.00	22.0
Bioreactor 1	1169.24	22.1
Bioreactor 1	1175.48	22.2
Bioreactor 1	1181.72	22.3
Bioreactor 1	1187.96	22.4
Bioreactor 1	1194.20	22.5
Bioreactor 1	1200.44	22.6
Bioreactor 1	1206.68	22.7
Bioreactor 1	1212.92	22.8
Bioreactor 1	1219.16	22.9
Bioreactor 1	1225.40	23.0
Bioreactor 1	1231.64	23.1
Bioreactor 1	1237.88	23.2
Bioreactor 1	1244.12	23.3
Bioreactor 1	1250.36	23.4
Bioreactor 1	1256.60	23.5
Bioreactor 1	1262.84	23.6
Bioreactor 1	1269.08	23.7
Bioreactor 1	1275.32	23.8
Bioreactor 1	1281.56	23.9
Bioreactor 1	1287.80	24.0
Bioreactor 1	1294.04	24.1
Bioreactor 1	1300.28	24.2
Bioreactor 1	1306.52	24.3
Bioreactor 1	1312.76	24.4
Bioreactor 1	1319.00	24.5
Bioreactor 1	1325.24	24.6
Bioreactor 1	1331.48	24.7
Bioreactor 1	1337.72	24.8
Bioreactor 1	1343.96	24.9
Bioreactor 1	1350.20	25.0
Bioreactor 1	1356.44	25.1
Bioreactor 1	1362.68	25.2
Bioreactor 1	1368.92	25.3
Bioreactor 1	1375.16	25.4
Bioreactor 1	1381.40	25.5
Bioreactor 1	1387.64	25.6
Bioreactor 1	1393.88	25.7
Bioreactor 1	1400.12	25.8
Bioreactor 1	1406.36	25.9
Bioreactor 1	1412.60	26.0
Bioreactor 1	1418.84	26.1
Bioreactor 1	1425.08	26.2
Bioreactor 1	1431.32	26.3
Bioreactor 1	1437.56	26.4
Bioreactor 1	1443.80	26.5
Bioreactor 1	1450.04	26.6
Bioreactor 1	1456.28	26.7
Bioreactor 1	1462.52	26.8
Bioreactor 1	1468.76	26.9
Bioreactor 1	1475.00	27.0
Bioreactor 1	1481.24	27.1
Bioreactor 1	1487.48	27.2
Bioreactor 1	1493.72	27.3
Bioreactor 1	1500.00	27.4
Bioreactor 1	1506.24	27.5
Bioreactor 1	1512.48	27.6
Bioreactor 1	1518.72	27.7
Bioreactor 1	1525.00	27.8
Bioreactor 1	1531.24	27.9
Bioreactor 1	1537.48	28.0
Bioreactor 1	1543.72	28.1
Bioreactor 1	1550.00	28.2
Bioreactor 1	1556.24	28.3
Bioreactor 1	1562.48	28.4
Bioreactor 1	1568.72	28.5
Bioreactor 1	1575.00	28.6
Bioreactor 1	1581.24	28.7
Bioreactor 1	1587.48	28.8
Bioreactor 1	1593.72	28.9
Bioreactor 1	1600.00	29.0
Bioreactor 1	1606.24	29.1
Bioreactor 1	1612.48	29.2
Bioreactor 1	1618.72	29.3
Bioreactor 1	1625.00	29.4
Bioreactor 1	1631.24	29.5
Bioreactor 1	1637.48	29.6
Bioreactor 1	1643.72	29.7
Bioreactor 1	1650.00	29.8
Bioreactor 1	1656.24	29.9
Bioreactor 1	1662.48	30.0
Bioreactor 1	1668.72	30.1
Bioreactor 1	1675.00	30.2
Bioreactor 1	1681.24	30.3
Bioreactor 1	1687.48	30.4
Bioreactor 1	1693.72	30.5
Bioreactor 1	1700.00	30.6
Bioreactor 1	1706.24	30.7
Bioreactor 1	1712.48	30.8
Bioreactor 1	1718.72	30.9
Bioreactor 1	1725.00	31.0
Bioreactor 1	1731.24	31.1
Bioreactor 1	1737.48	31.2
Bioreactor 1	1743.72	31.3
Bioreactor 1	1750.00	31.4
Bioreactor 1	1756.24	31.5
Bioreactor 1	1762.48	31.6
Bioreactor 1	1768.72	31.7
Bioreactor 1	1775.00	31.8
Bioreactor 1	1781.24	31.9
Bioreactor 1	1787.48	32.0
Bioreactor 1	1793.72	32.1
Bioreactor 1	1800.00	32.2
Bioreactor 1	1806.24	32.3
Bioreactor 1	1812.48	32.4
Bioreactor 1	1818.72	32.5
Bioreactor 1	1825.00	32.6
Bioreactor 1	1831.24	32.7
Bioreactor 1	1837.48	32.8
Bioreactor 1	1843.72	32.9
Bioreactor 1	1850.00	33.0
Bioreactor 1	1856.24	33.1
Bioreactor 1	1862.48	33.2
Bioreactor 1	1868.72	33.3
Bioreactor 1	1875.00	33.4
Bioreactor 1	1881.24	33.5
Bioreactor 1	1887.48	33.6
Bioreactor 1	1893.72	33.7
Bioreactor 1	1900.00	33.8
Bioreactor 1	1906.24	33.9
Bioreactor 1	1912.48	34.0
Bioreactor 1	1918.72	34.1
Bioreactor 1	1925.00	34.2
Bioreactor 1	1931.24	34.3
Bioreactor 1	1937.48	34.4
Bioreactor 1	1943.72	34.5
Bioreactor 1	1950.00	34.6
Bioreactor 1	1956.24	34.7
Bioreactor 1	1962.48	34.8
Bioreactor 1	1968.72	34.9
Bioreactor 1	1975.00	35.0
Bioreactor 1	1981.24	35.1
Bioreactor 1	1987.48	35.2
Bioreactor 1	1993.72	35.3
Bioreactor 1	2000.00	35.4
Bioreactor 1	2006.24	35.5
Bioreactor 1	2012.48	35.6
Bioreactor 1	2018.72	35.7
Bioreactor 1	2025.00	35.8
Bioreactor 1	2031.24	35.9
Bioreactor 1	2037.48	36.0
Bioreactor 1	2043.72	36.1
Bioreactor 1	2050.00	36.2
Bioreactor 1	2056.24	36.3
Bioreactor 1	2062.48	36.4
Bioreactor 1	2068.72	36.5
Bioreactor 1	2075.00	36.6
Bioreactor 1	2081.24	36.7
Bioreactor 1	2087.48	36.8
Bioreactor 1	2093.72	36.9
Bioreactor 1	2100.00	37.0
Bioreactor 1	2106.24	37.1
Bioreactor 1	2112.48	37.2
Bioreactor 1	2118.72	37.3
Bioreactor 1	2125.00	37.4
Bioreactor 1	2131.24	37.5
Bioreactor 1	2137.48	37.6
Bioreactor 1	2143.72	37.7
Bioreactor 1	2150.00	37.8
Bioreactor 1	2156.24	37.9
Bioreactor 1	2162.48	38.0
Bioreactor 1	2168.72	38.1
Bioreactor 1	2175.00	38.2
Bioreactor 1	2181.24	38.3
Bioreactor 1	2187.48	38.4
Bioreactor 1	2193.72	38.5
Bioreactor 1	2200.00	38.6
Bioreactor 1	2206.24	38.7
Bioreactor 1	2212.48	38.8
Bioreactor 1	2218.72	38.9
Bioreactor 1	2225.00	39.0
Bioreactor 1	2231.24	39.1
Bioreactor 1	2237.48	39.2
Bioreactor 1	2	

# DOE vs MVDA for Ambr<sup>®</sup> Experiments

## MODDE<sup>®</sup> DOE

Media Component Screening

Cell Line and Clone Screening

Bioreactor Condition Optimization

Feed Strategy Optimization

Bioreactor Characterization

Scale-up/down Model Development

VS

## SIMCA<sup>®</sup> MVDA

Bioreactor Condition Optimization

Feed Strategy Optimization

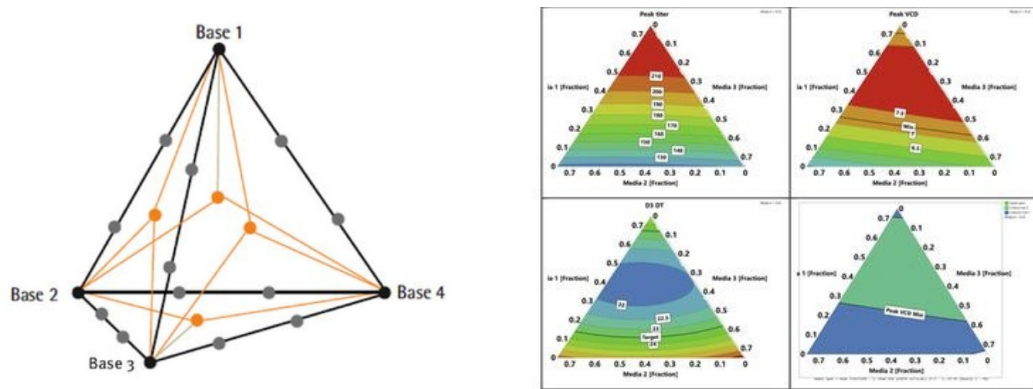
Spectroscopy Calibration Modeling

Scale-up/down Model Verification



# Ambr<sup>®</sup> DOE Examples

## Media Component Screening



### Challenge:

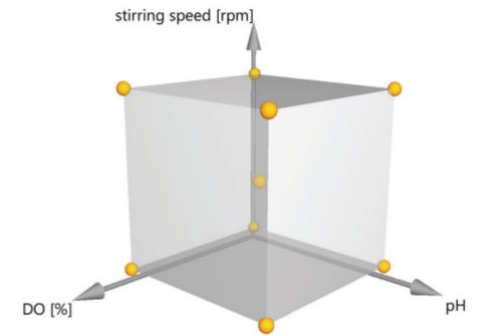
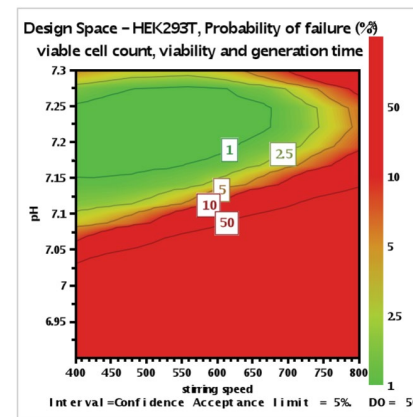
- Not all commercially available media supports every cell line, therefore in some cases additional supplements must be added to improve cell performance
- The classical approach to media component screening is time consuming and labor-intensive

### Data Driven Solution:

- Use DOE principals to systematically screen media components for productivity effects
- Recognize interactions and define them early in the process, setting optimum levels accordingly

How to Optimize Cell Culture Media to Speed Biopharma Development, Sartorius (2018)

## Bioreactor Optimization and Characterization



### Challenge:

- A significant challenge for developing viral vector gene therapies is that is well characterized and can be scaled-up

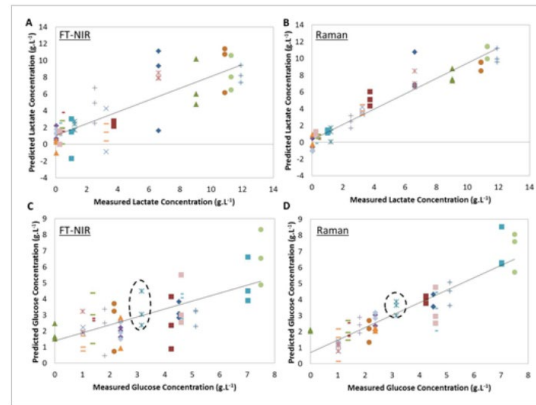
### Data Driven Solution

- Use DOE principals to identify which process parameters impact product quality and yield
- Justify and adjust manufacturing operating ranges - control strategy - and acceptance criteria

Optimization of HEK293t Suspension Cultivation with DOE-approach in the Ambr 15 Micro Bioreactor, Bollman, Riethmuller, Johansson, Tappe, R&D Regenerative Medicine, Sartorius (2019)

# Ambr<sup>®</sup> MVDA Examples

## Spectroscopy Calibration Modeling



### Challenge

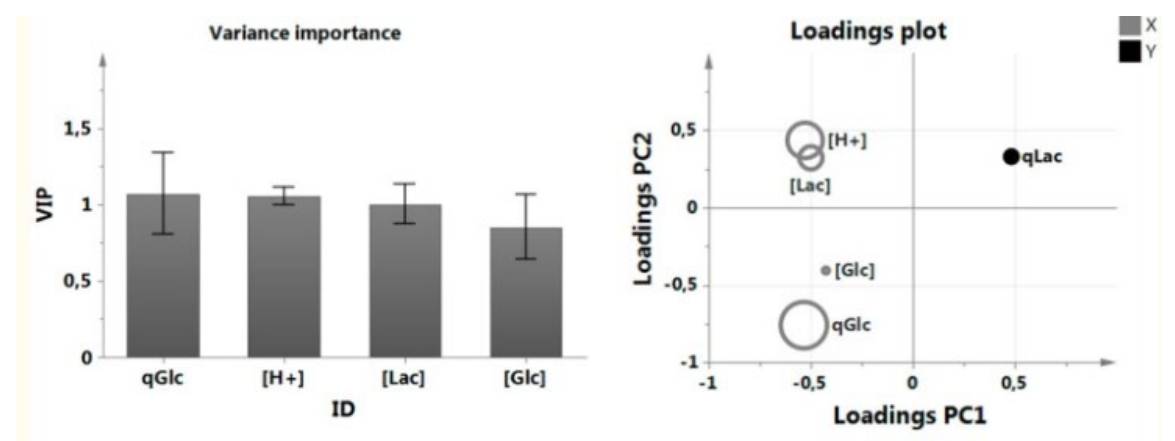
- PAT tools can (i.e. spectral devices, advanced sensors and analyzers) generate large amounts of complex data presenting a challenge when it comes to interpreting accuracy

### Data Driven Solution:

- MVDA is a fast and flexible spectral calibration tool that can handle multiple types of spectral data (NIR, IR, Raman, Fluorescence, Mass-Spec)
- It provides accurate prediction models for analyte concentrations can be used to optimize monitoring and control of the culture

Comparison of Spectroscopy Technologies for Improved Monitoring of Cell Culture Processes in Miniature Bioreactors, Rowland-Jones, van den Berg, Racher, Martin, Jaques, Biotechnology Process Volume 33, Issue 2, GSK (2017)

## Feed-Strategy Optimization



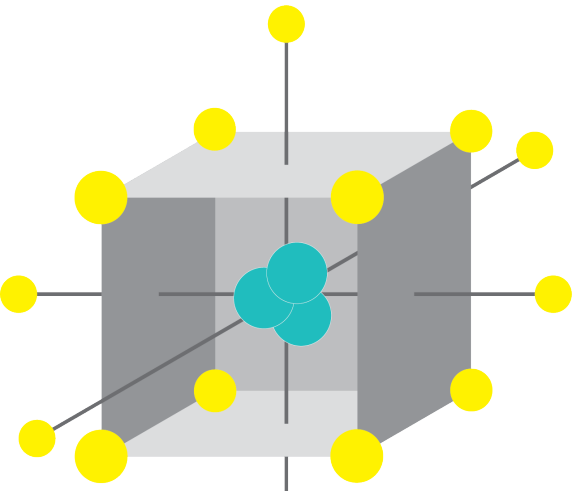
### Challenge

- A significant challenge in cell-culture cultivation is predicting the nutritional requirements of the culture so that an appropriate feeding strategy can be implemented
- Data Driven Solution
- Use MVDA to identify and rank any potential inhibitors or promoters of cell-culture productivity in order to optimize feed strategy

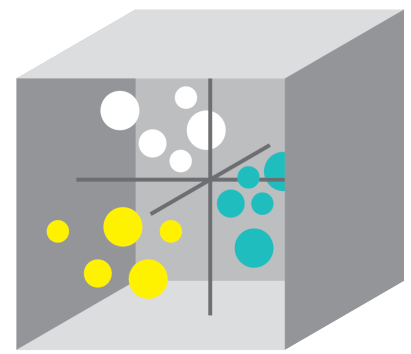
Metabolic Control in Mammalian Fed-Batch Cell Cultures for Reduced Lactic Acid Accumulation and Improved Process Robustness, Kanokovsky, Clemens, Müller, Bechmann, Berger, Schlatter, Herwig, Boehringer Ingelheim, Vienna Institute of Technology, Bioengineering Basel (2016)



# Combining the Best of Both Worlds for Ambr<sup>®</sup> Scale-down Modeling



DOE  
Design of Experiments



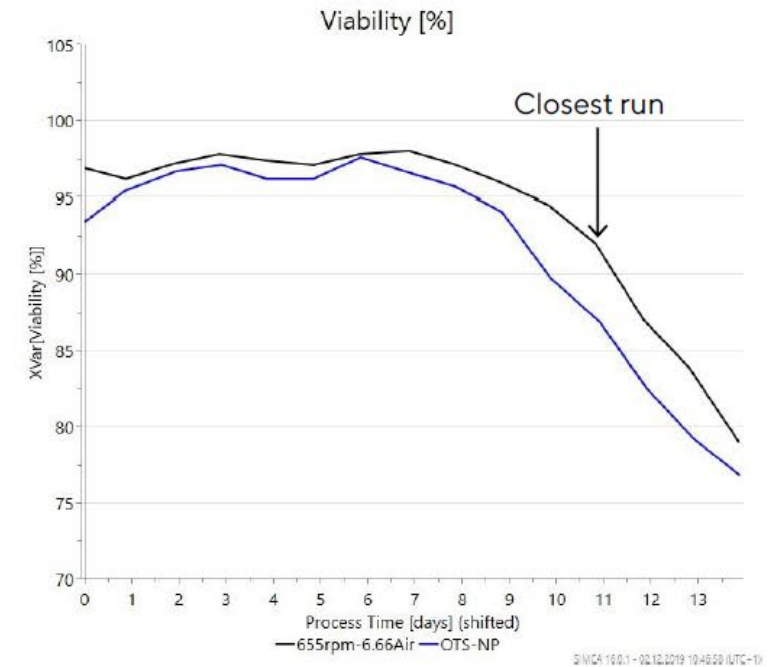
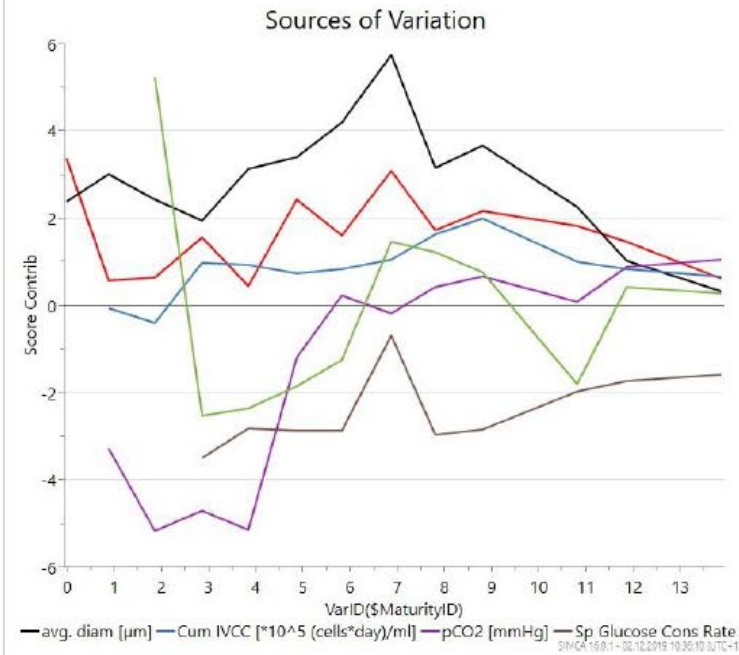
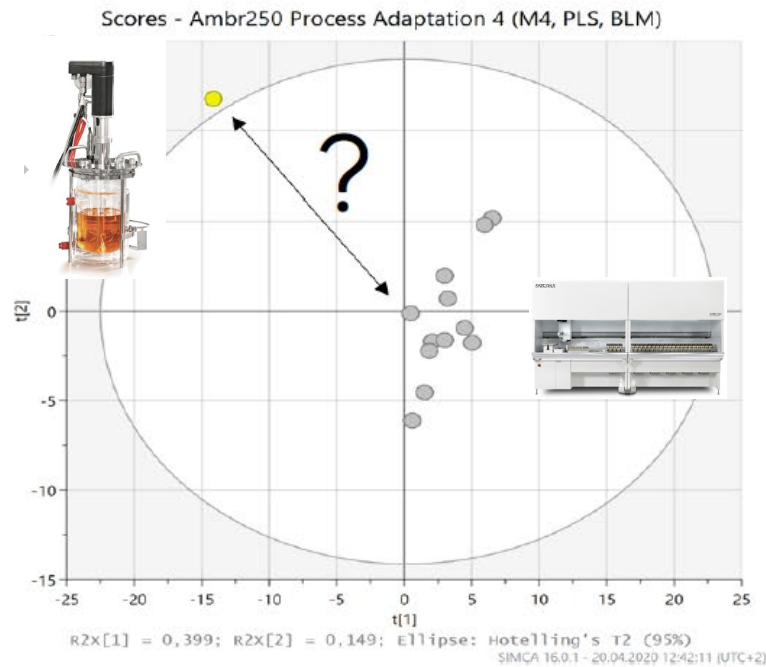
MVDA  
Multivariate Data Analysis



Combine multi-objective optimization criteria (DOE) with quantitative ranking (MVDA)

Squeeze all the value from the data you have

# DOE and MVDA for Ambr<sup>®</sup> Scale-down Modeling

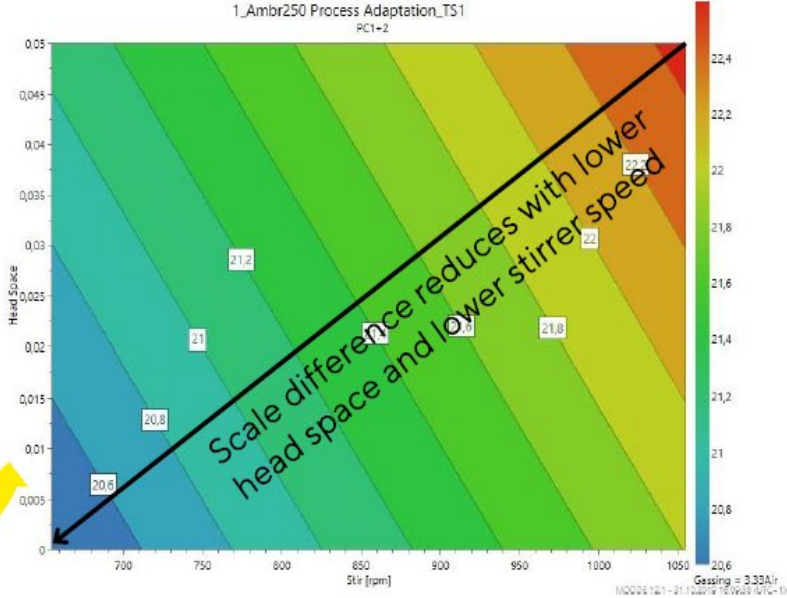
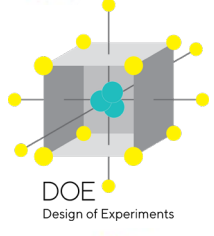
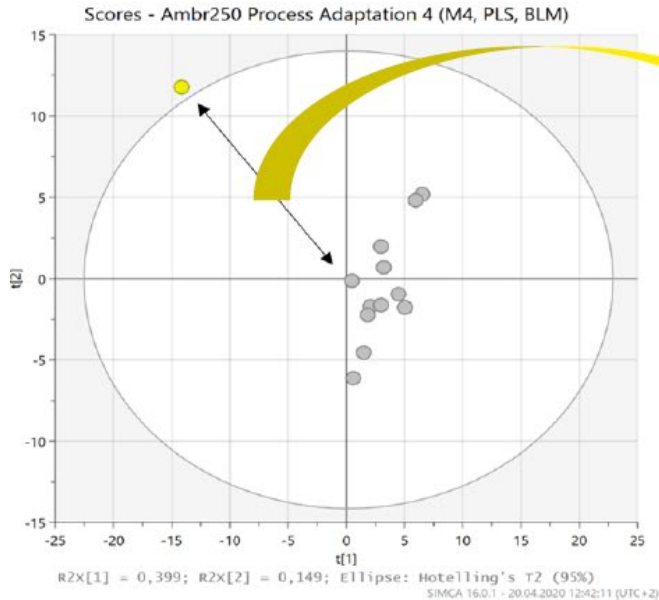


MVDA Used to Evaluate Major Sources of Variability Between Ambr<sup>®</sup> and 5L Reference Scale

DOE Establishment of Scale-down Bioprocess Models for Ambr Using Clustered Multivariate Analysis, Timo Schmidberger, Thomas Krieg, Sartorius (2020)

# DOE and MVDA for Ambr<sup>®</sup> Scale-down Modeling

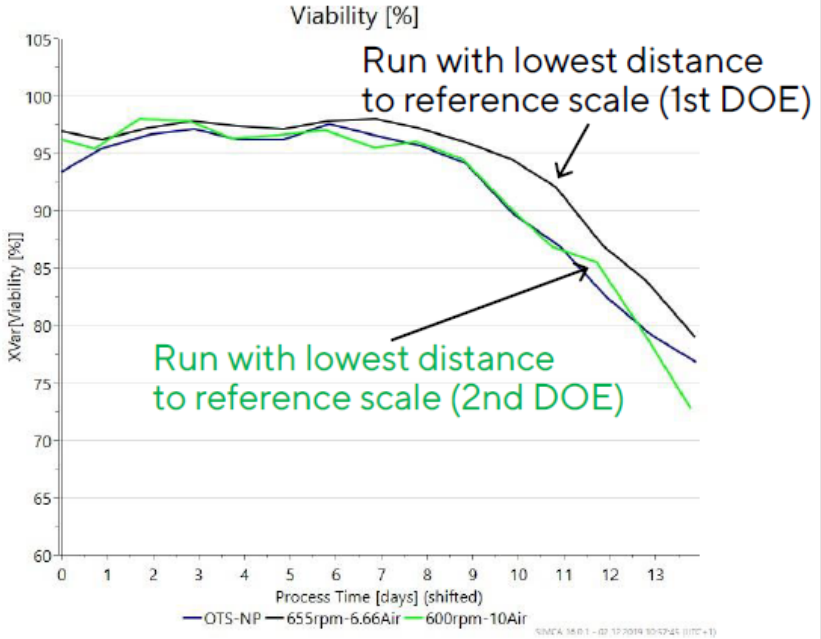
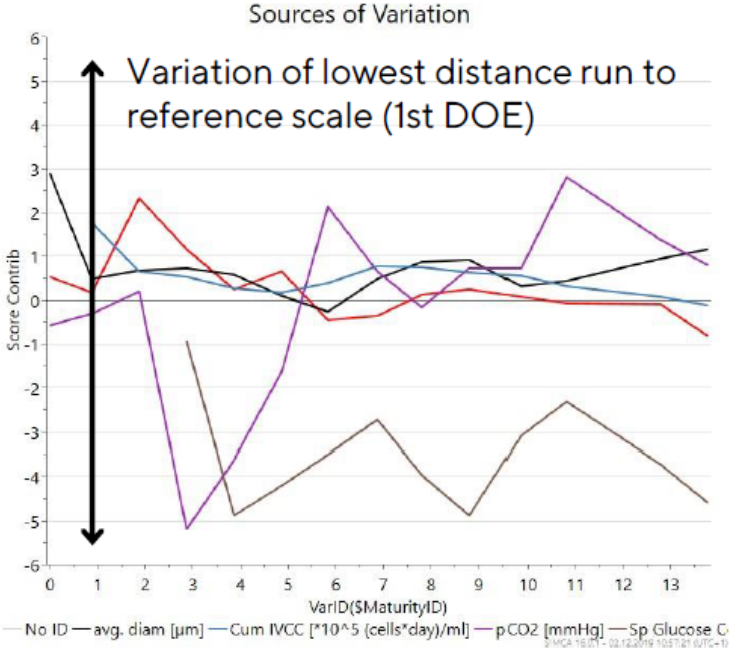
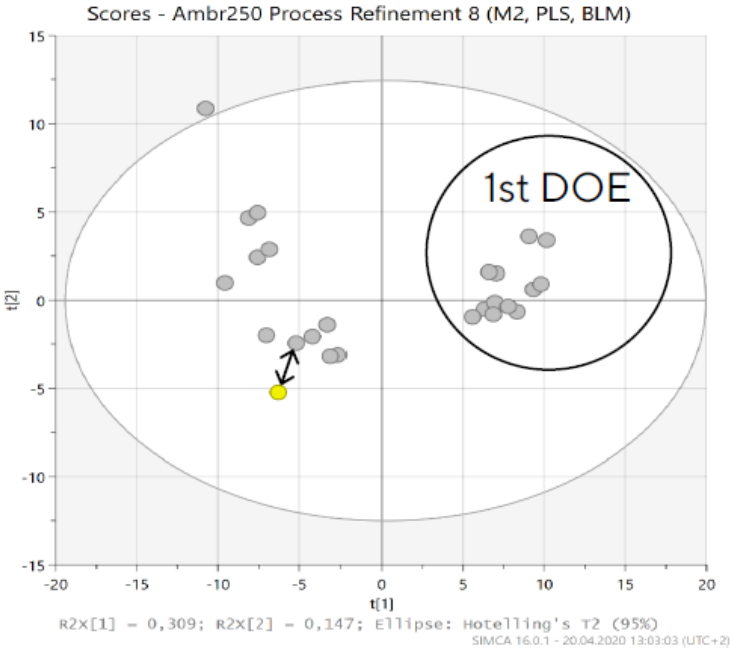
- The goal in scale-down modeling is to decrease the distance between the scaled-down runs and the target scale
- To do this we can use MODDE<sup>®</sup> DOE to evaluate the length  $\longleftrightarrow$  and give suggestions as to how to change operating parameters to reduce scale difference



DOE Establishment of Scale-down Bioprocess Models for Ambr Using Clustered Multivariate Analysis, Timo Schmidberger, Thomas Krieg, Sartorius (2020)



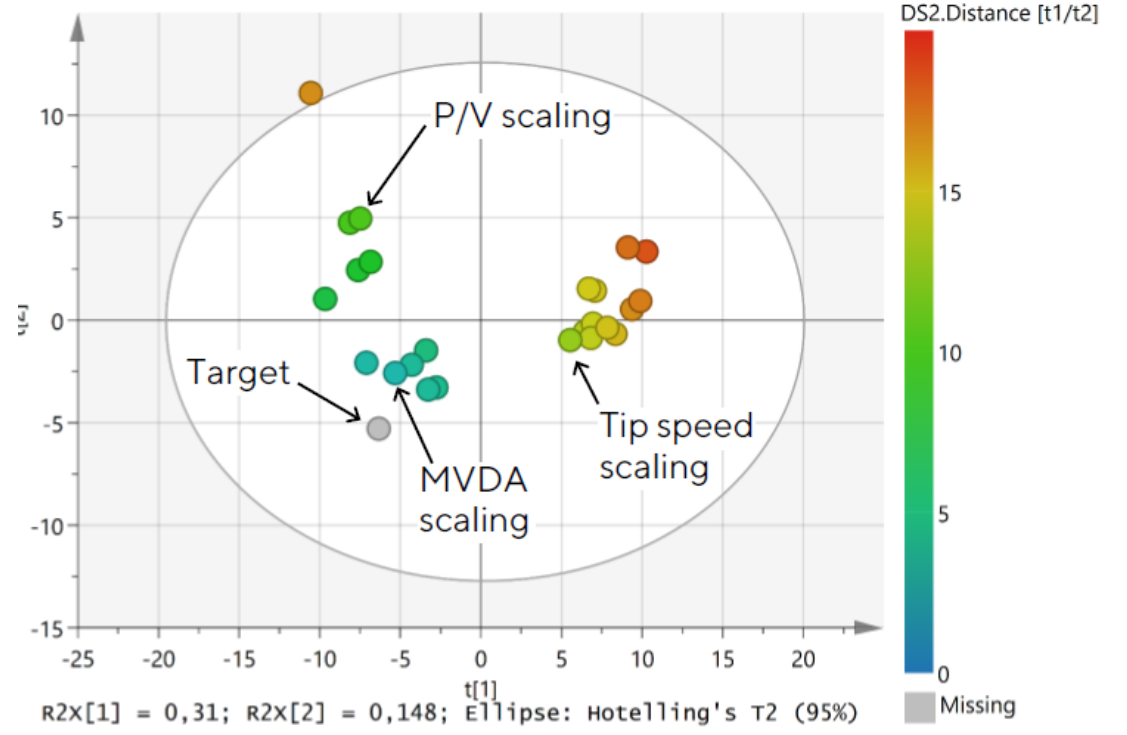
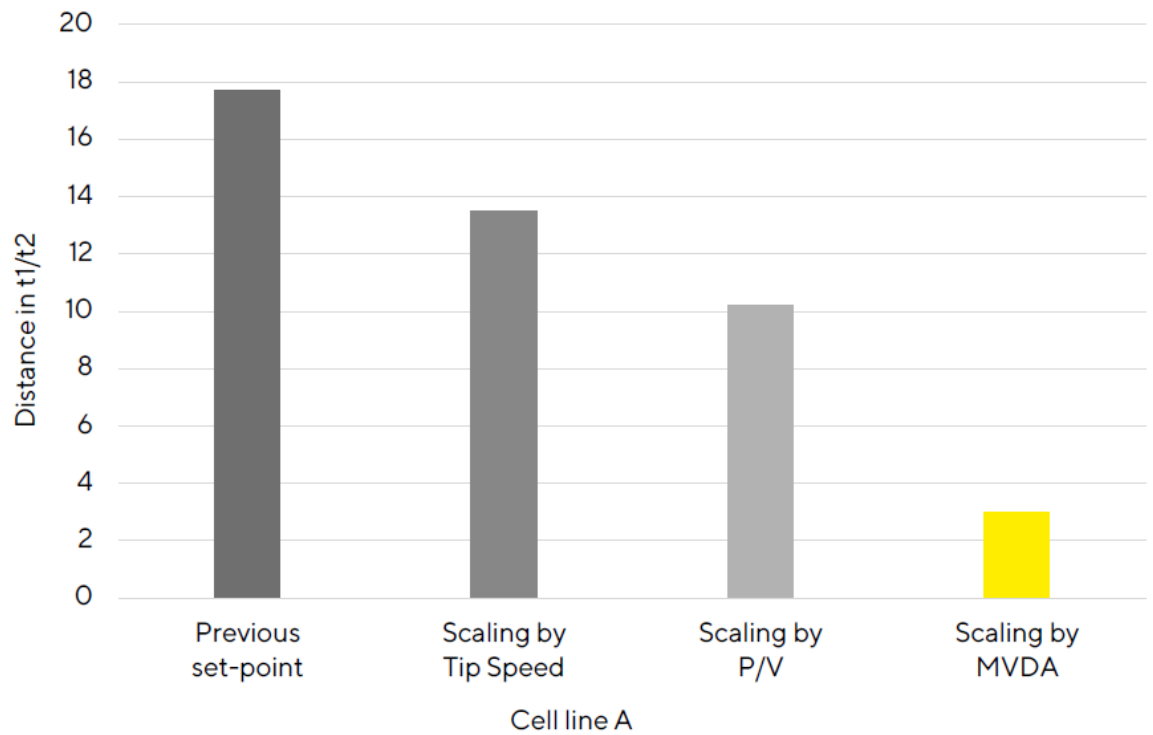
# DOE and MVDA for Ambr<sup>®</sup> Scale-down Modeling



DOE Establishment of Scale-down Bioprocess Models for Ambr Using Clustered Multivariate Analysis, Timo Schmidberger, Thomas Krieg, Sartorius (2020)



# DOE and MVDA for Ambr<sup>®</sup> Scale-down Modeling Method Conformation



DOE Establishment of Scale-down Bioprocess Models for Ambr Using Clustered Multivariate Analysis, Timo Schmidberger, Thomas Krieg, Sartorius (2020)

# DOE and QbD are Key for CDMO Success

“The most successful CDMOs, therefore, have identified strategies for completing process development projects efficiently and effectively while incorporating DOE and QbD approaches that provide increased process understanding and lead to optimal processes.”

Greg Flyte, GSK, CMO Alliance and Program Management





Visit the CDMO Landing page

[www.landing.umetrics.com/en/cdmo](http://www.landing.umetrics.com/en/cdmo)



Read the latest blogs and watch the recorded webinars



Try the ROI calculator



Book a strategy session

# Visit Sartorius at IFPAC 2021



Join Presentation on Hybrid Modeling for Deeper Bioprocess Insight by Catalina Moreno on March 3

IFPAC 2021 Digital Event



Join Presentation on Bringing Bioprocess Digital Twins to Life by Tiffany McLeod and Chris McCready on March 1

IFPAC 2021 Digital Event



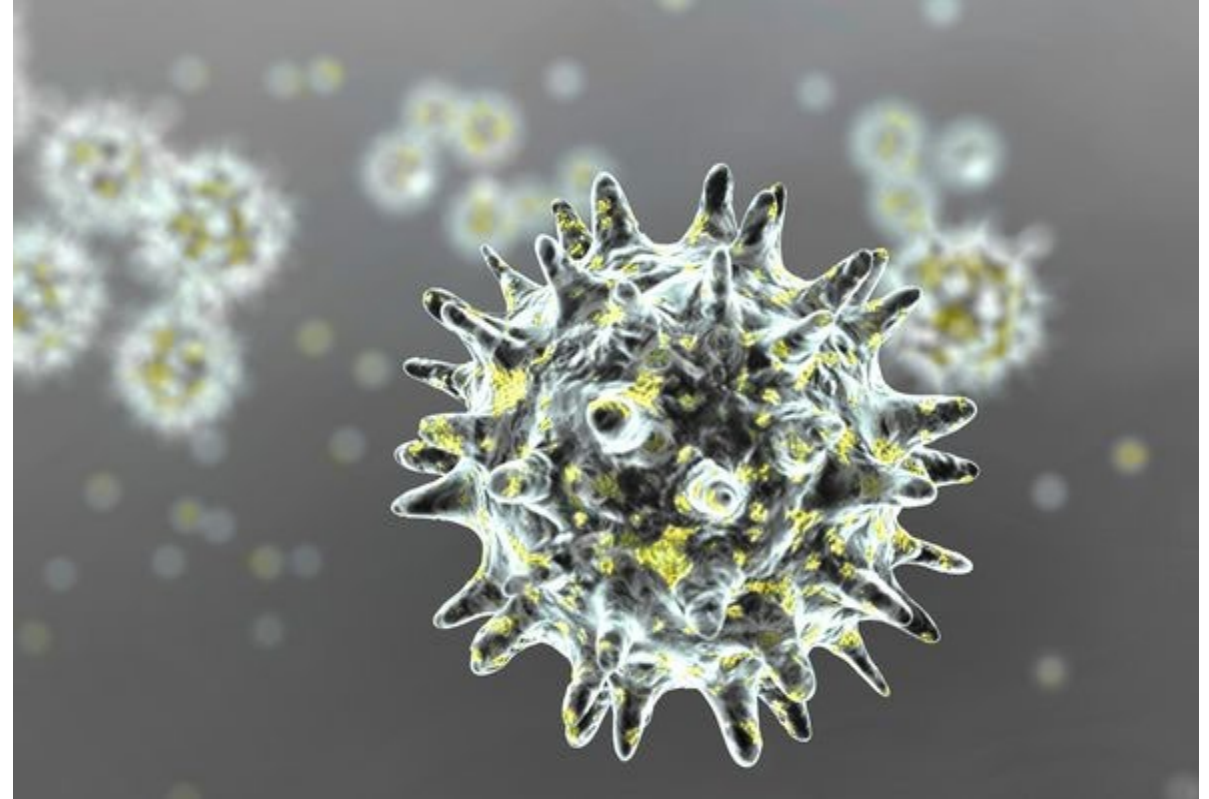
# How CDMOs can Digitalize their Cell and Gene Therapy Processes

Join our next webinar in the CDMO Webinar Series

March 24<sup>th</sup> 4-5 PM CET | 10-11 AM EST

More Info Can Be Found on the CDMO landing page

[www.landing.umetrics.com/en/cdmo](http://www.landing.umetrics.com/en/cdmo)



Thank you!

Tiffany McLeod  
Kevin McHugh

**SARTORIUS**