

Simplifying Progress

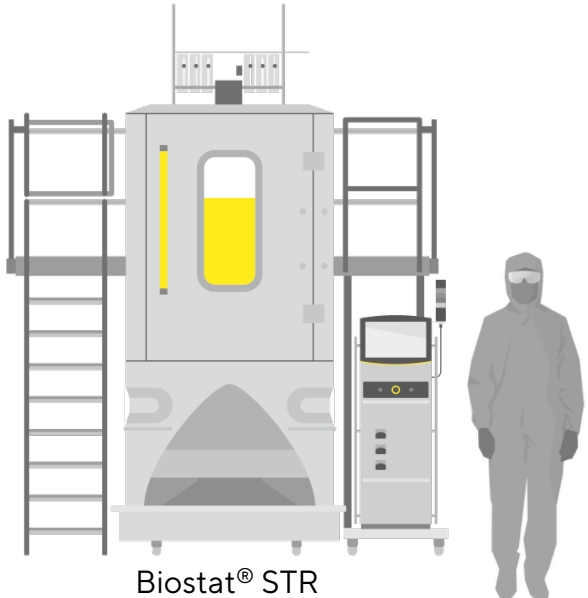
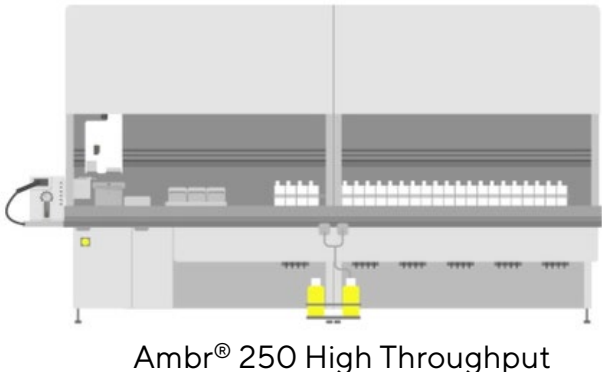
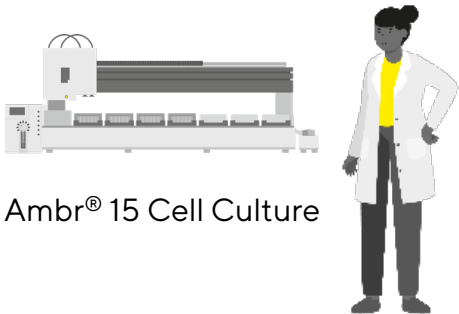


## Introducing BioPAT<sup>®</sup> Spectro: Unlocking the Potential of Raman Spectroscopy

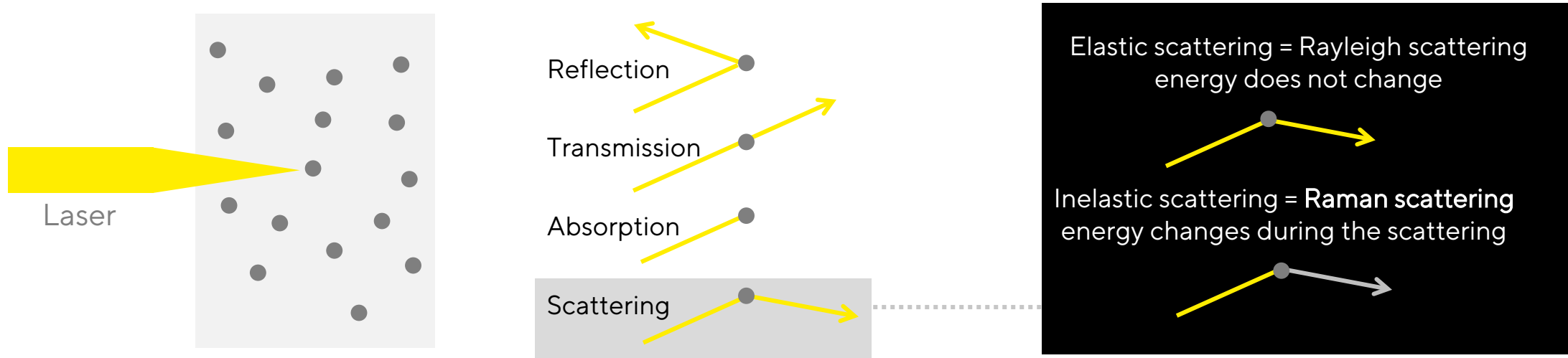
Veronica Brivio - PAT Product Manager, Dan Kopec - PAT Specialist for North America | September 2022

**SARTORIUS**

# BioPAT<sup>®</sup> Spectro Introduces Raman Spectroscopy Into Ambr<sup>®</sup> and Biostat STR<sup>®</sup> to Enable Quality by Design (QbD)



# Raman Is a Laser Scattering Technique With High Molecular Specificity



A laser is projected into the bioreactor

- Non-invasive
- In-line

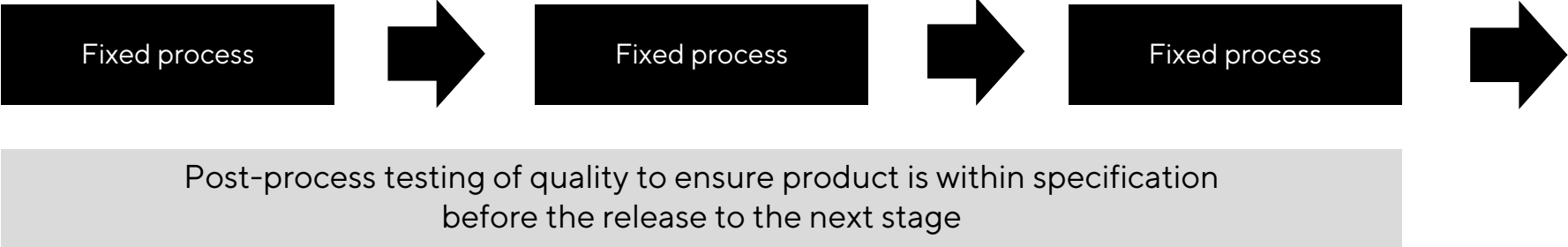
The laser interacts with molecules in a non-destructive fashion

Inelastic scattering is measured as a Raman spectrum

- Gives “molecular fingerprint”, enabling qualitative and quantitative analysis of several analytes, such as glucose, lactate, titer etc
- Weak signal requires longer acquisition time

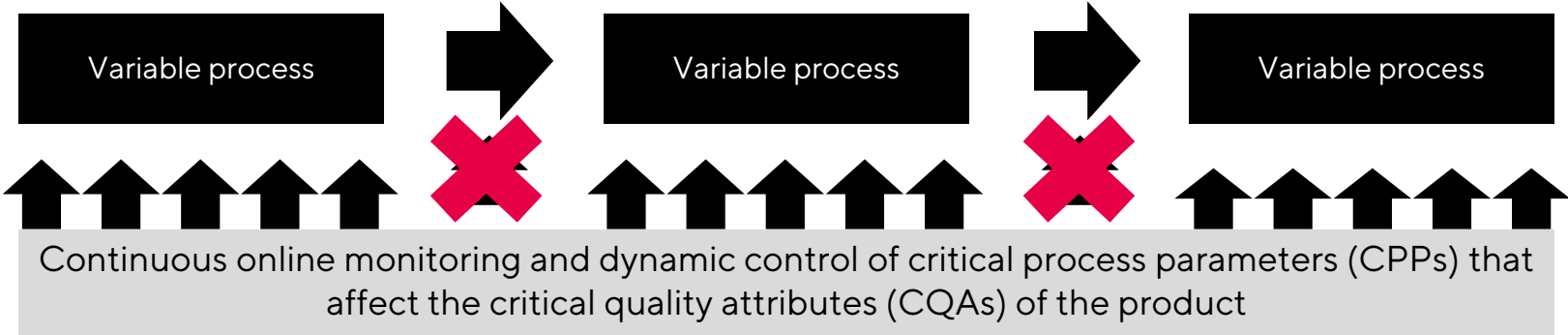
# Quality by Design Enables Consistent and High Product Quality

## Traditional process – Post-process quality testing



Variability in product quality and quantity

## PAT process – Quality by Design (QbD)



Consistency in product quality and quantity

Real-time release

# Raman Spectroscopy Is Complex and Not Well Integrated With Single-Use



Expectation



Reality



## Process development

Full DOE and model building is too laborious and expensive  
missing automation and high throughput integrations



## Data management

Data from different sources require manual alignment  
requires expert data scientist and is error prone



## SU manufacturing

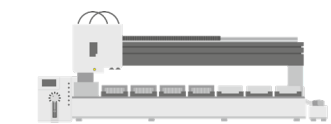
No comprehensive single-use integration  
immersion probes suffer from stray light in transparent bags

Limited scalability due to low-robustness models and lack of scalable hardware

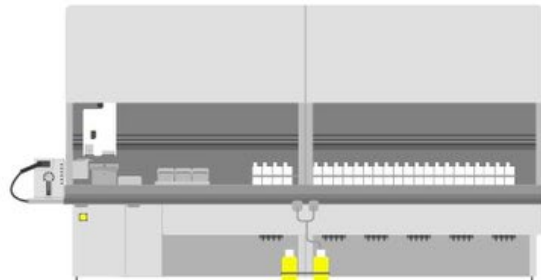
# BioPAT® Spectro Was Designed to Meet Three Key Requirements

- Enable Raman spectroscopy in high throughput process development
- Facilitate and improve the model building and data management process
- Full single-use integration and scalability for commercial manufacturing

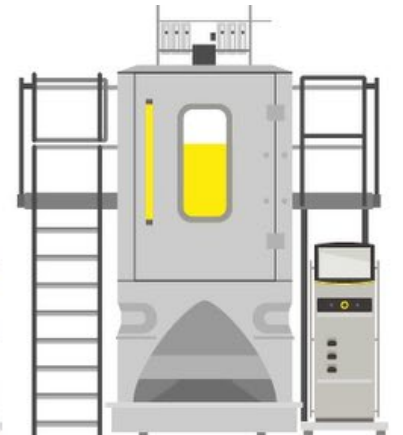
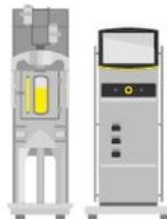
'[BioPAT® Spectro is] the next killer app for Raman as it enables the technology to expand beyond manufacturing and into PD.'



Ambr® 15 Cell Culture



Ambr® 250 High Throughput

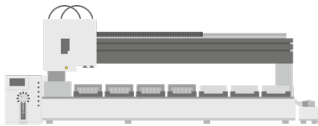


Biostat® STR

Aspen Alert, March 2020

# Raman Spectroscopy Use Cases

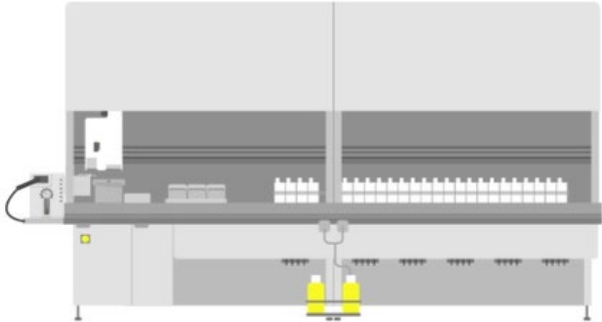
Model building



Ambr® 15 Cell Culture

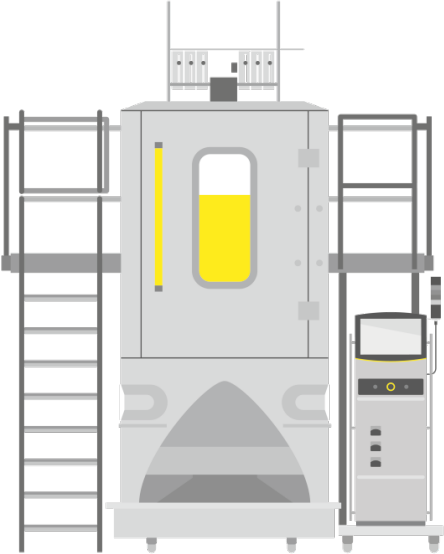


Process development



Ambr® 250 High Throughput

Single-use manufacturing



Biostat® STR

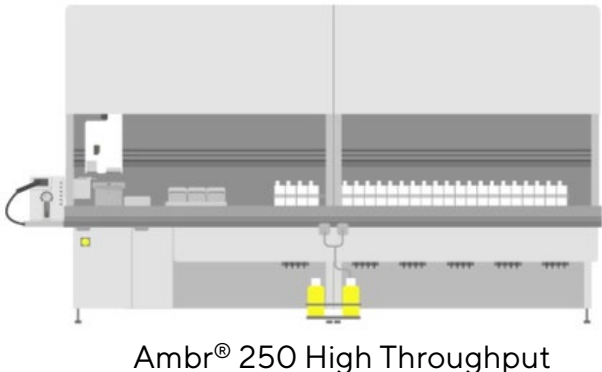
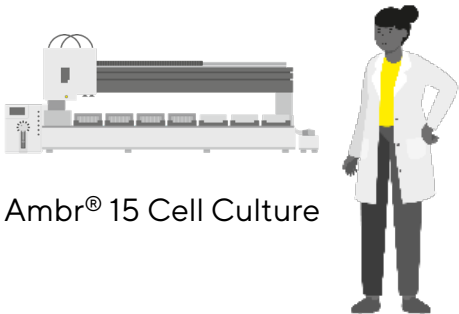


# Raman Spectroscopy Use Cases

Model building

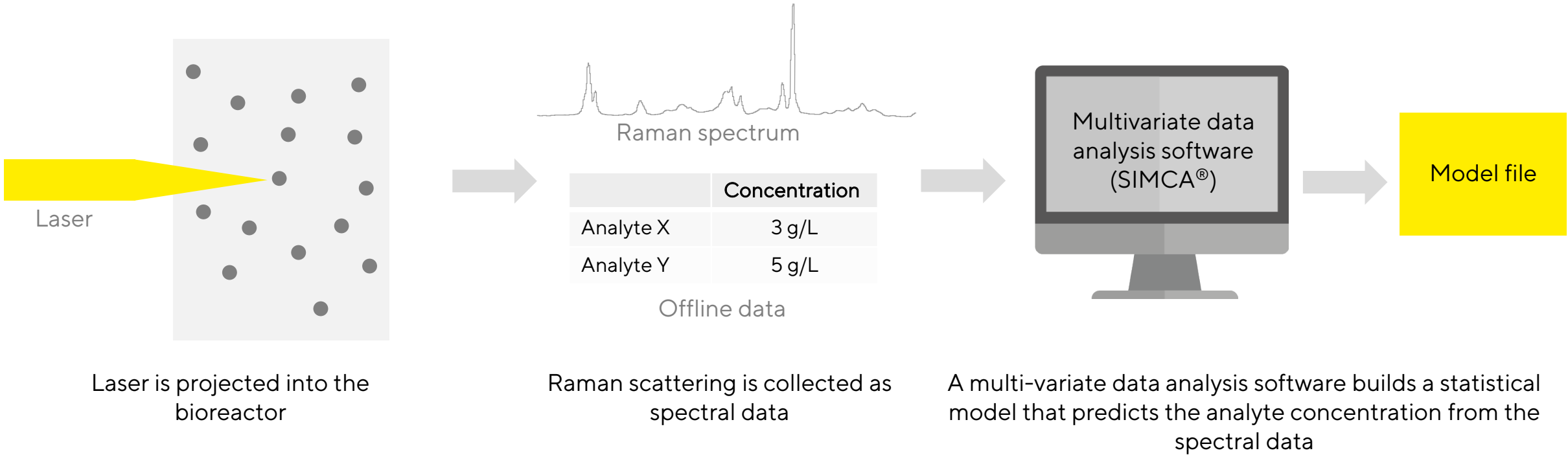
Process development

Single-use manufacturing

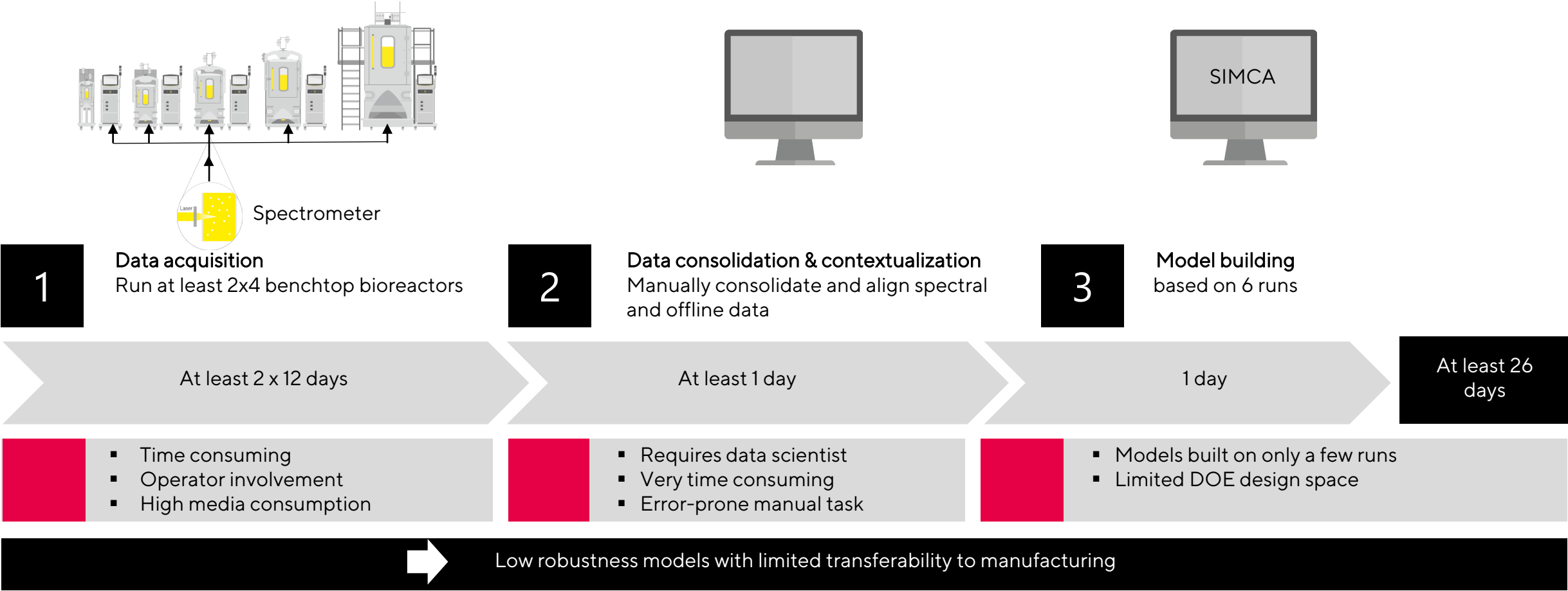




# Quantitative Raman Spectroscopy Requires Model Building



# Conventional Model Building Is Laborious, Expensive and Time-Consuming



# BioPAT<sup>®</sup> Spectro is integrated in Ambr<sup>®</sup> 15 and Ambr<sup>®</sup> 250 High Throughput

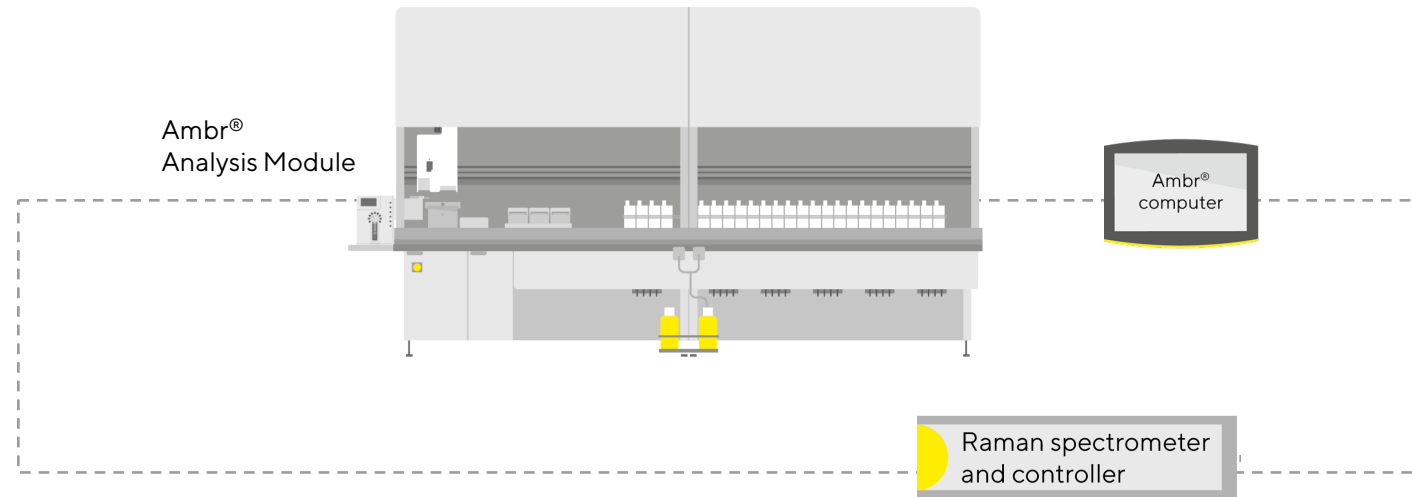


Connection of a Endress+Hauser Raman spectrometer probe



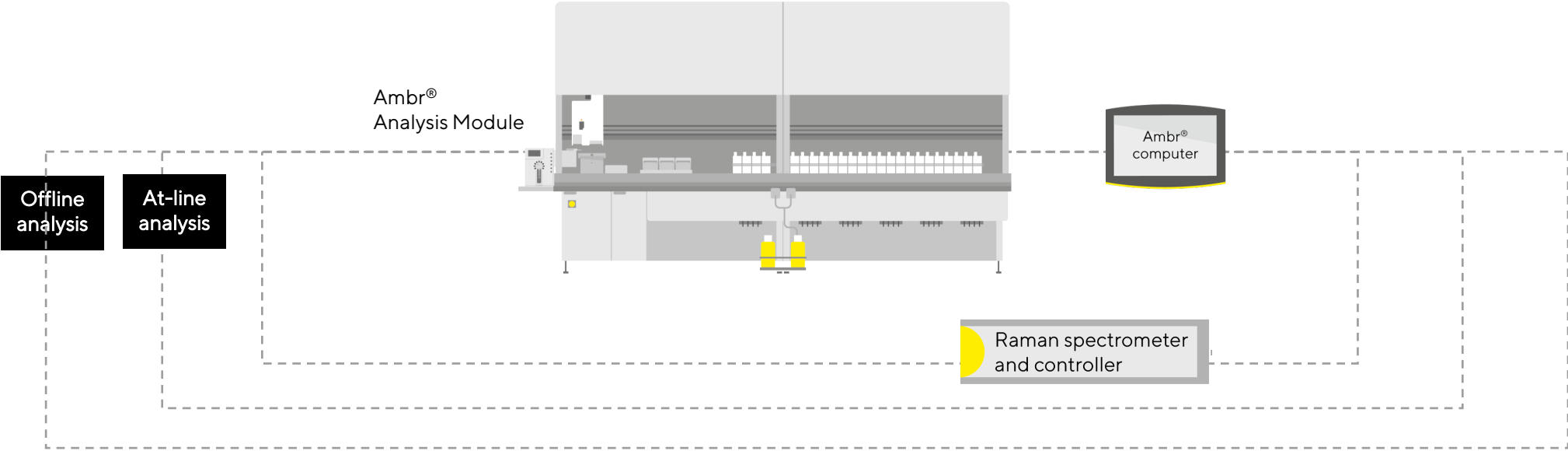
Connection of a Tornado Raman spectrometer probe

# Data Acquisition and Consolidation Is Fully Automated in Ambr<sup>®</sup>



3<sup>rd</sup> party Raman spectrometers (Kaiser Optical Systems and Tornado Spectral Systems) are fully integrated into the Ambr<sup>®</sup> and can be controlled via the Ambr<sup>®</sup> software. Spectral data is collected in the Ambr<sup>®</sup> software.

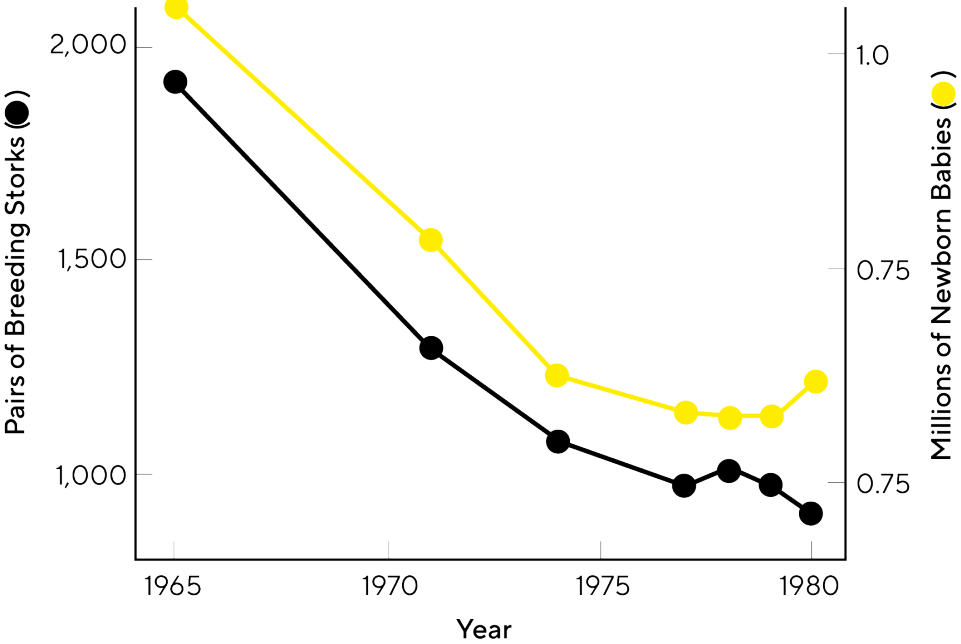
# Data Acquisition and Consolidation Is Fully Automated in Ambr®



The data from integrated analyzers is collected and automatically aligned with the spectral data.  
Offline data can be added manually during the run.



# BioPAT<sup>®</sup> Spectro in Ambr<sup>®</sup> Is Designed for Building Causal Models

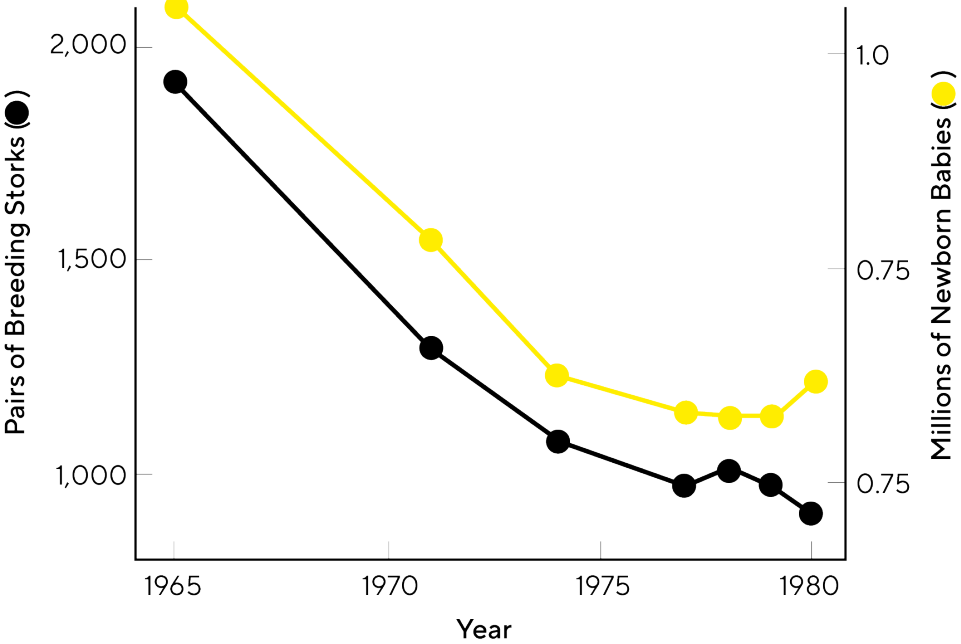


SIES, H.A new parameter for sex education. Nature 223. 495 (1988)

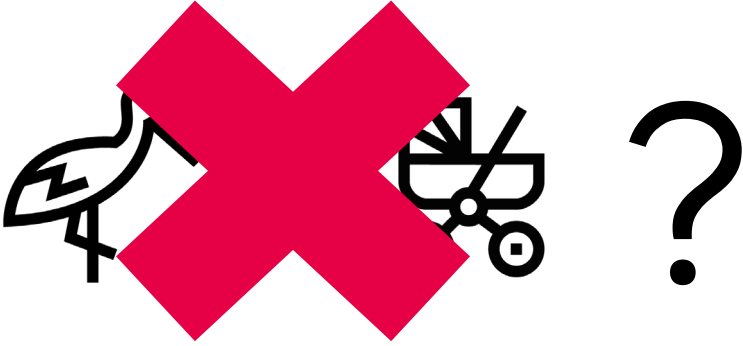


Image: © John Lund/Getty Images

# BioPAT<sup>®</sup> Spectro in Ambr<sup>®</sup> Is Designed for Building Causal Models



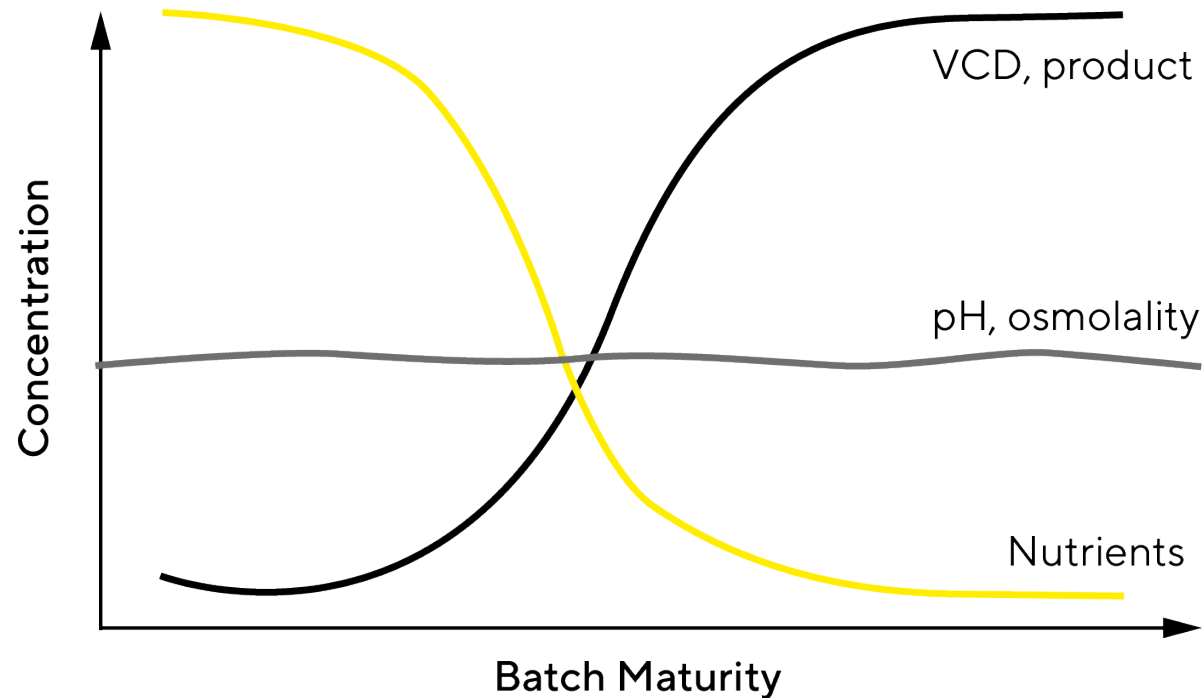
SIES, H.A new parameter for sex education. Nature 223. 495 (1988)



Correlation does not imply causation!



# BioPAT<sup>®</sup> Spectro in Ambr<sup>®</sup> Is Designed for Building Causal Models



## Random correlations are a risk in model building

- Analyte concentration changes follow similar patterns
- statistical models cannot distinguish causal from non-causal correlations
- Risk mitigation: Use techniques to break correlations

# BioPAT® Spectro in Ambr® Is Designed for Building Causal Models

## Experimental plan

24 vessels in Ambr® 250 or Ambr® 15

18 vessel DOE breaks correlations and increases model robustness

6 golden batches

Inoculation density

Glucose set point / profile

pH

Dissolved oxygen

## Automated spiking

Measure (unspiked) sample

Robot takes a new sample and adds stock solution from a well plate and mixes

Measure spiked sample

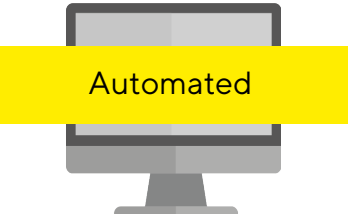
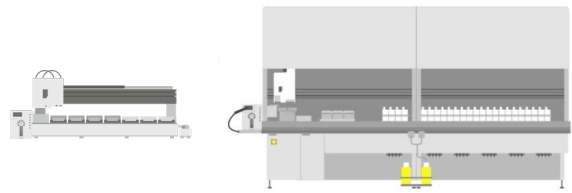
## Design of Experiment (DOE)

- Different process trajectories & analyte trends
- Trains model for predicting even during process deviations
- Breaks correlations & increases model robustness

## Analyte spiking

- Introduces a step change in concentration of an analyte
- Breaks correlations with non-spiked analytes
- Increases the concentration range of the model

# BioPAT<sup>®</sup> Spectro Is the Perfect Tool for Spectroscopy Model Building



1

## Data acquisition

Automated data acquisition & consolidation Run 12, 24 or 48 bioreactors in parallel

2

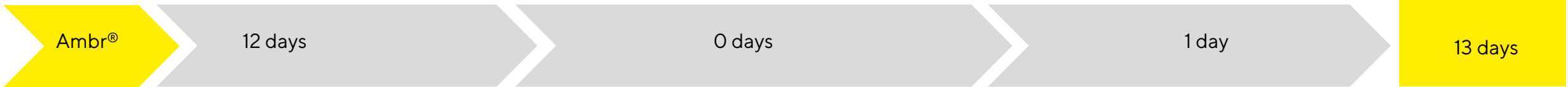
## Data consolidation & contextualization

Manually consolidate and align spectral and offline data

3

## Model building

Based on at least 10 runs



- Save 80% of operator time
- Save 95% of media costs
- Automated spiking breaks correlations

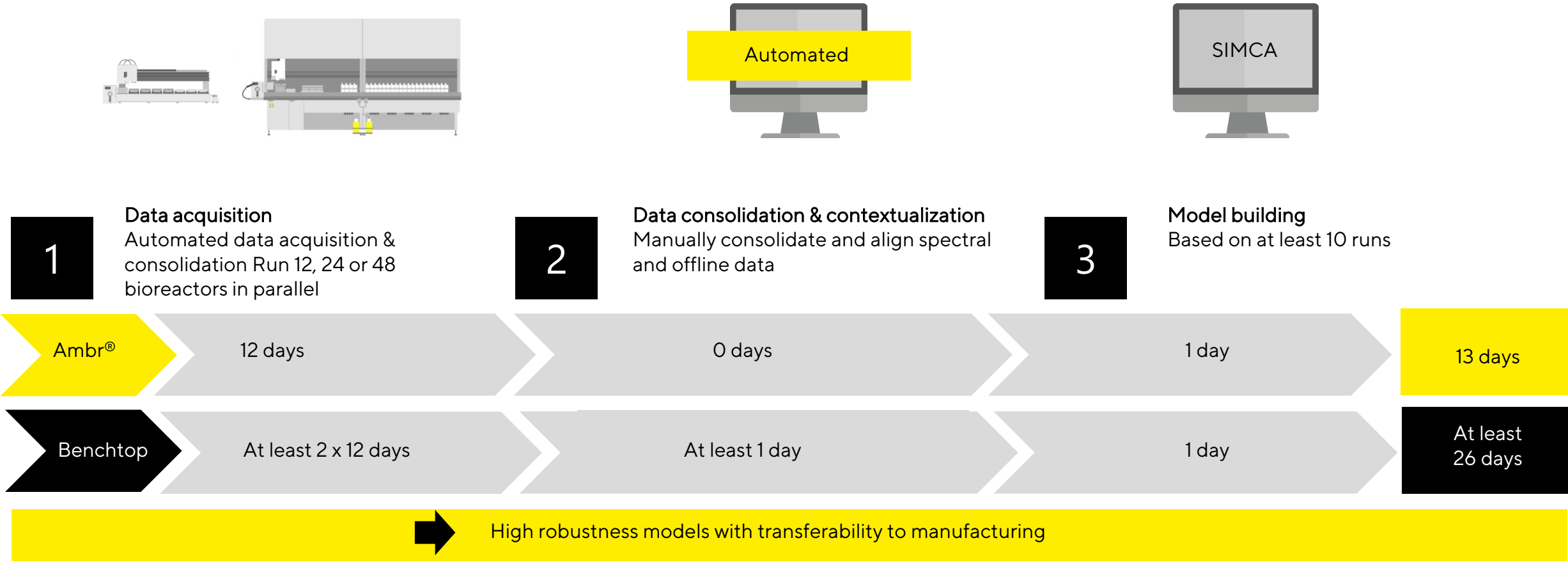
- Fully automated alignment of process, spectral and offline data
- Ready for important into SIMCA<sup>®</sup>

- Models built on many runs
- Large DOE design space
- All process data used

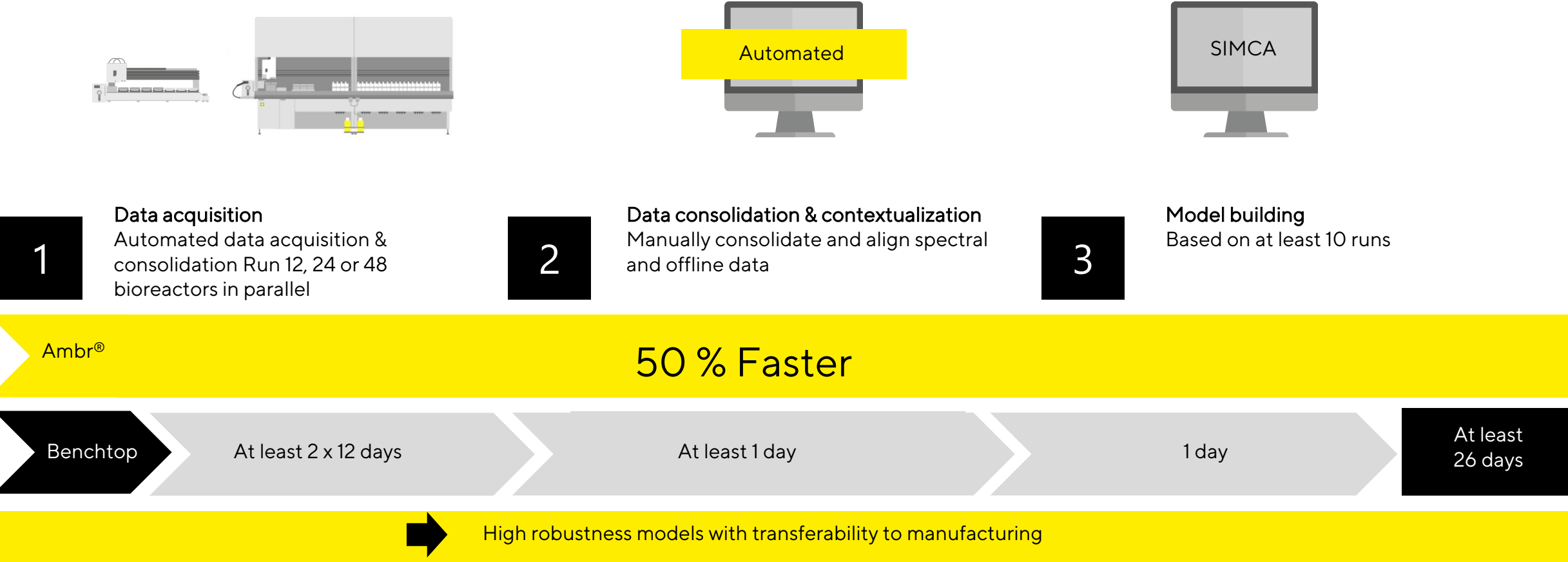


High robustness models with transferability to manufacturing

# BioPAT<sup>®</sup> Spectro Is the Perfect Tool for Spectroscopy Model Building



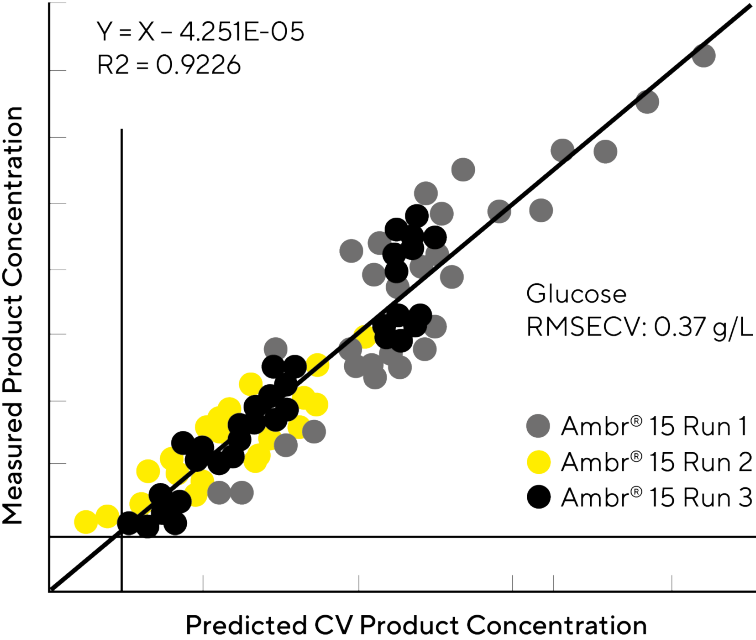
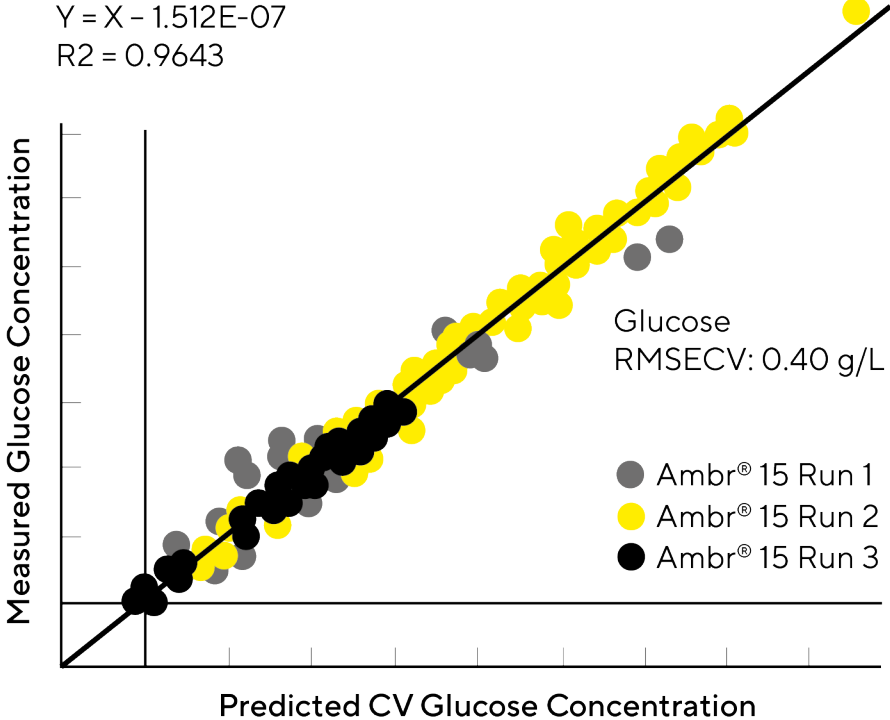
# BioPAT<sup>®</sup> Spectro Is the Perfect Tool for Spectroscopy Model Building



# Model Building In Ambr<sup>®</sup> 15 Cell Culture

## Predictive Models

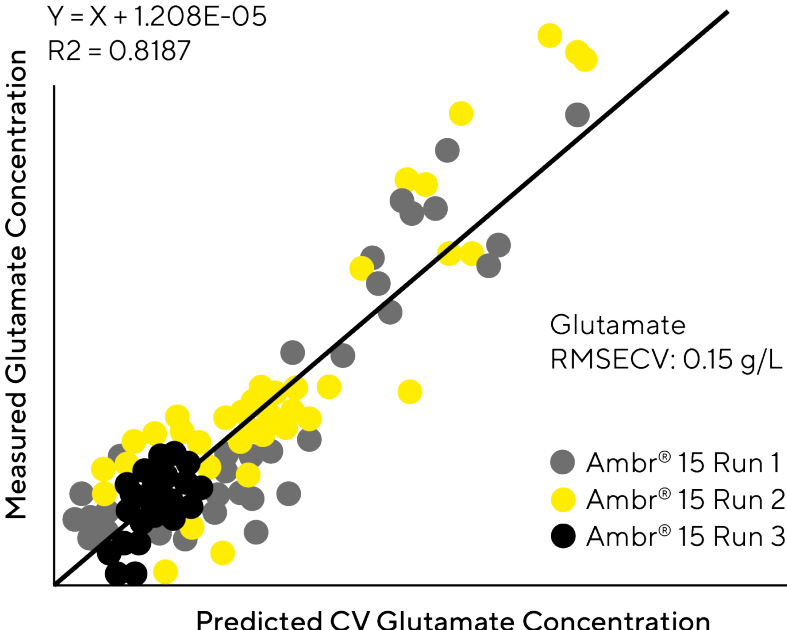
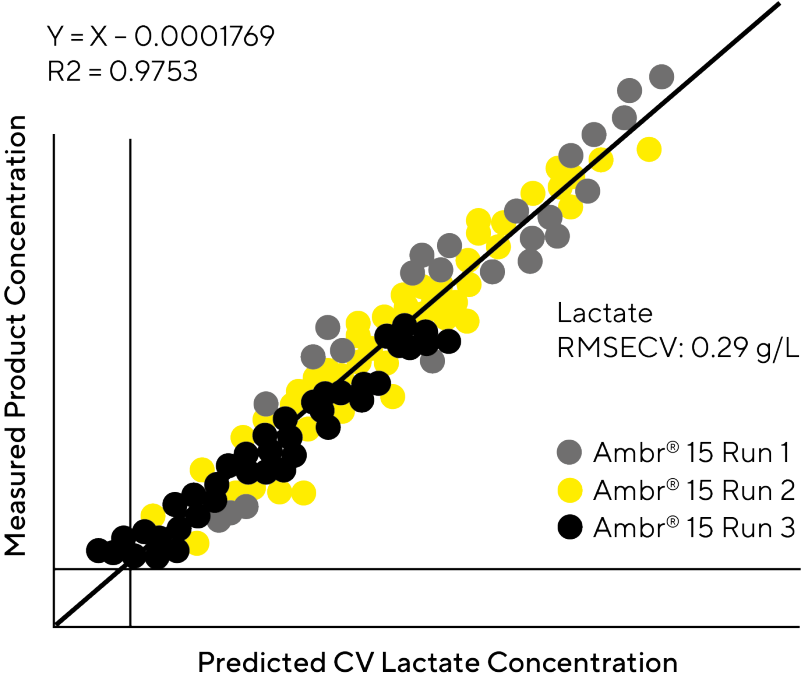
$Y = X - 1.512E-07$   
 $R^2 = 0.9643$



Reproduced with kind permission from Ruth Rowland-Jones, GSK Stevenage. Performed with a prototype spectroscopy integration and a Tornado Raman spectrometer.

# Model Building In Ambr<sup>®</sup> 15 Cell Culture

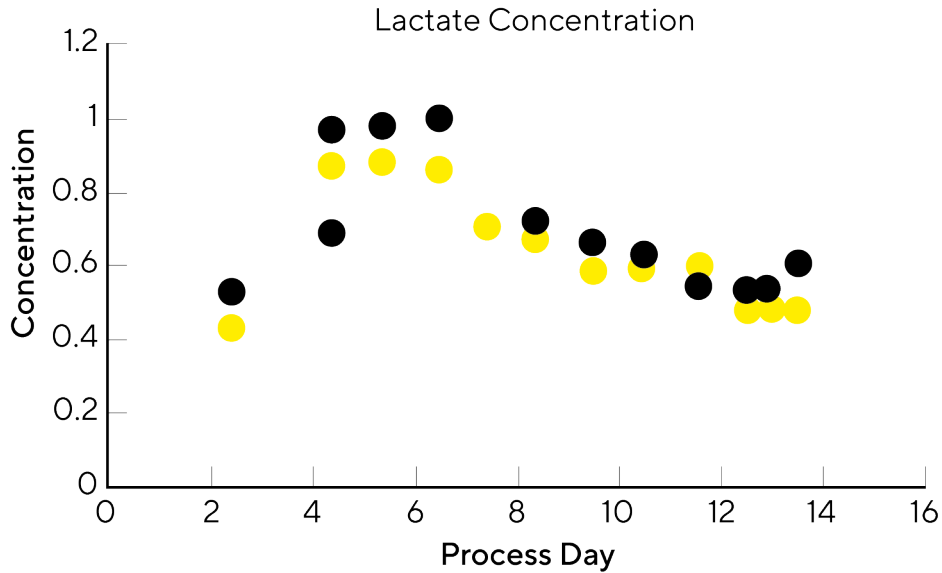
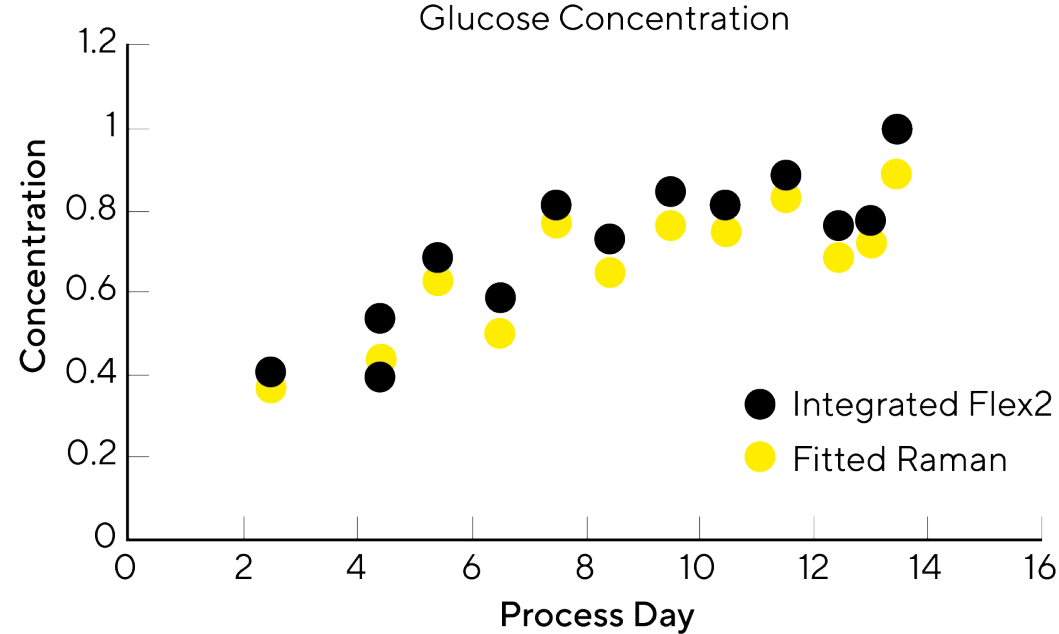
## Predictive Models



Reproduced with kind permission from Ruth Rowland-Jones, GSK Stevenage. Performed with a prototype spectroscopy integration and a Tornado Raman spectrometer.

# Model Building In Ambr<sup>®</sup> 15 Cell Culture

## Results - Single Bioreactor Time Course Comparison



All concentrations normalized to highest Flex2 Value

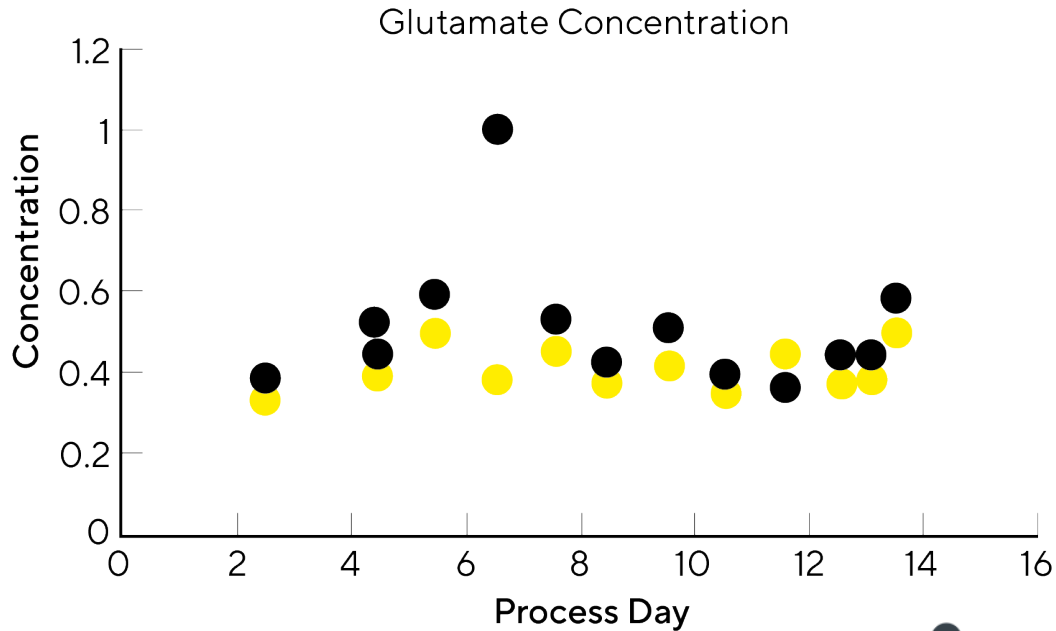
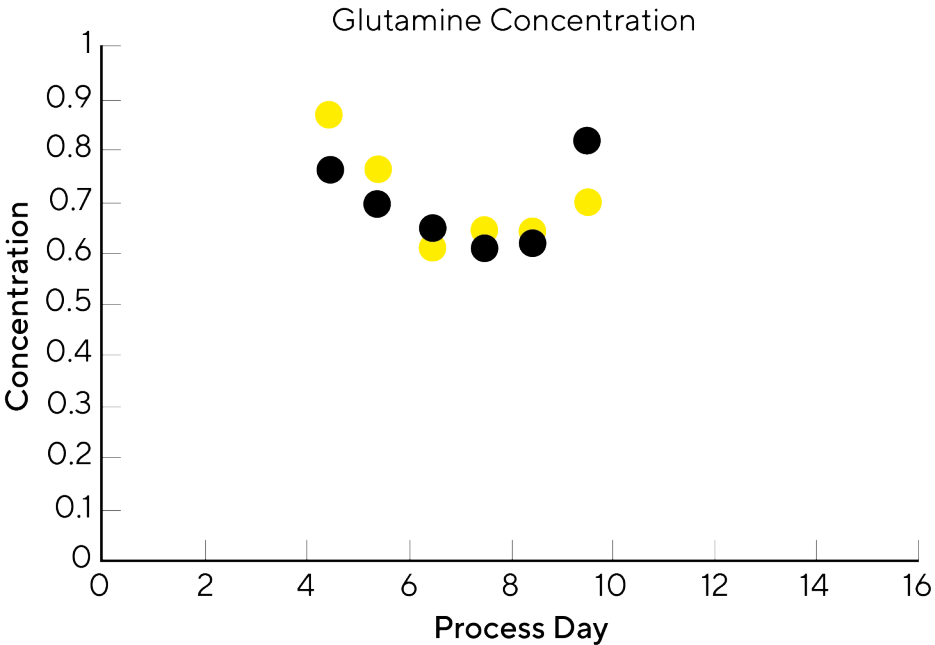


Reproduced with kind permission of Merck MSD, Mike Nelson presented this at AICHE 2019, performed with a prototype spectroscopy integration and a Tornado Raman spectrometer



# Model Building In Ambr<sup>®</sup> 15 Cell Culture

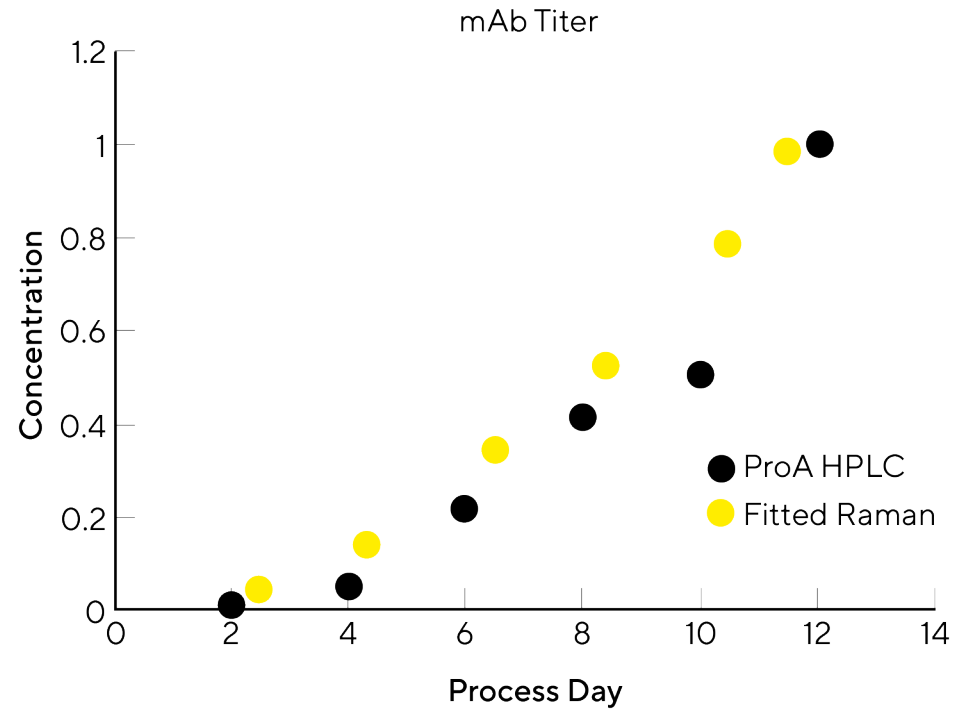
Results - Single Bioreactor Time Course Comparison



Reproduced with kind permission of Merck MSD, Mike Nelson presented this at AICHE 2019, performed with a prototype spectroscopy integration and a Tornado Raman spectrometer

# Model Building In Ambr<sup>®</sup> 15 Cell Culture

- Picked one specific bioreactor to show time course data
- Raman model tracks well with reference data
- Could conceivably control process using only fitted Raman model



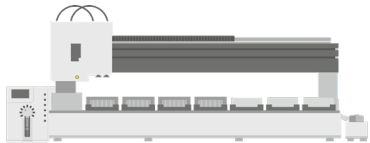
Reproduced with kind permission of Merck MSD, Mike Nelson presented this at AICHE 2019, performed with a prototype spectroscopy integration and a Tornado Raman spectrometer

# Raman Spectroscopy Use Cases

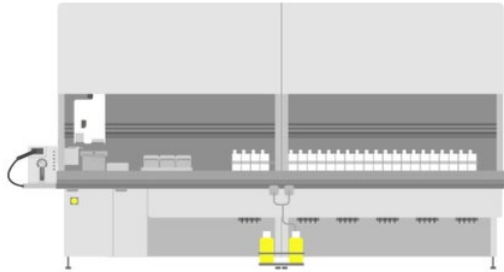
Model building

Process development

Single-use manufacturing

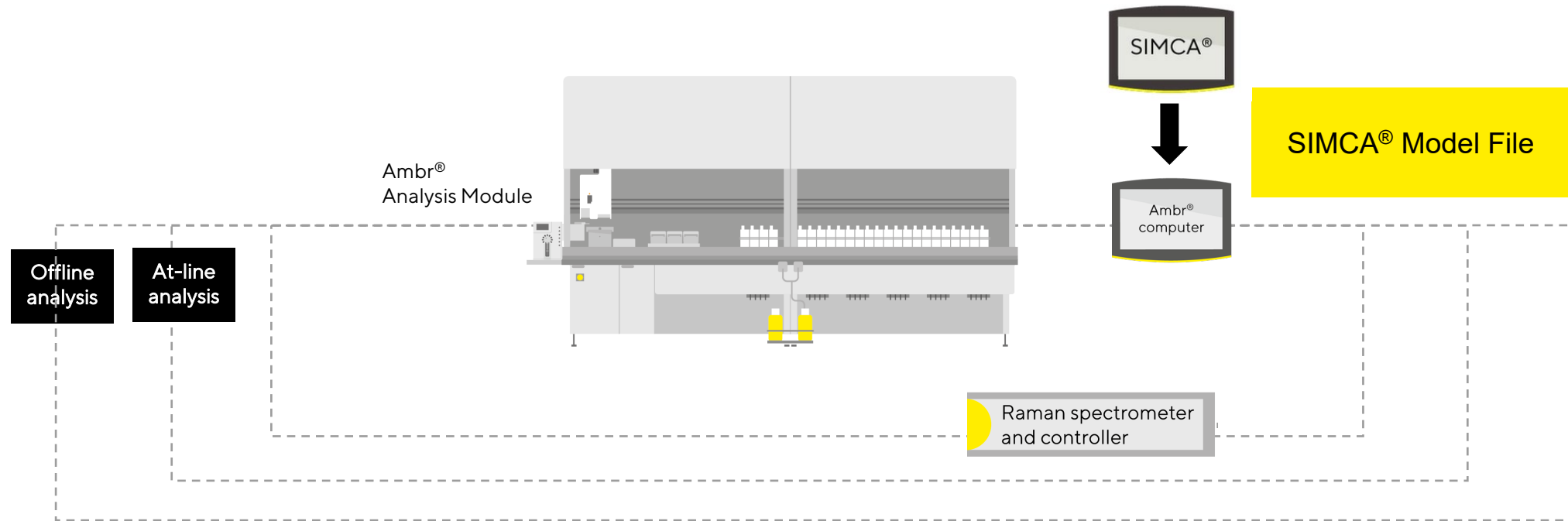


Ambr® 15 Cell Culture



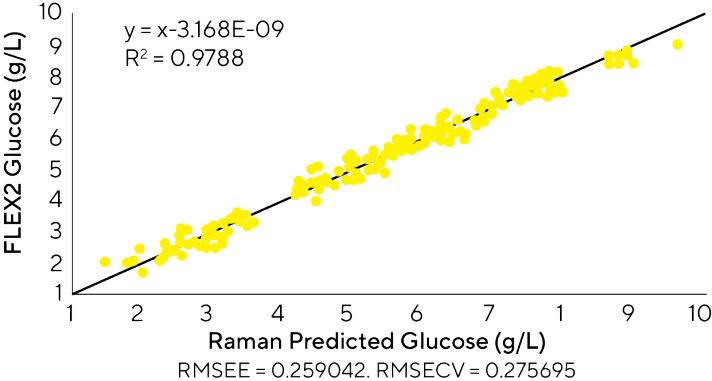
Ambr® 250 High Throughput

# Raman Spectroscopy in High-Throughput Process Development

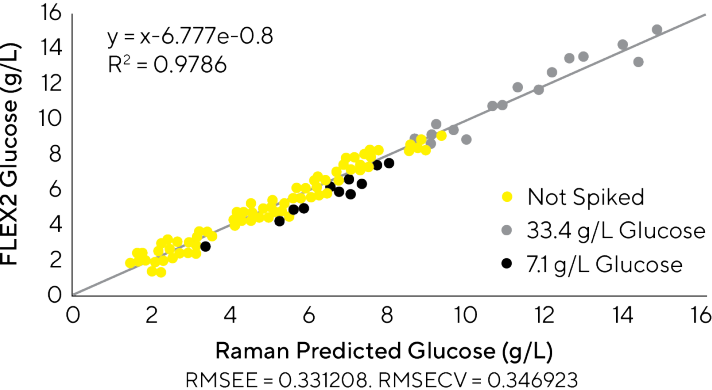


The Ambr<sup>®</sup> software is able to read SIMCA<sup>®</sup> model files and can predict analyte concentrations from the spectral data for process monitoring and control.

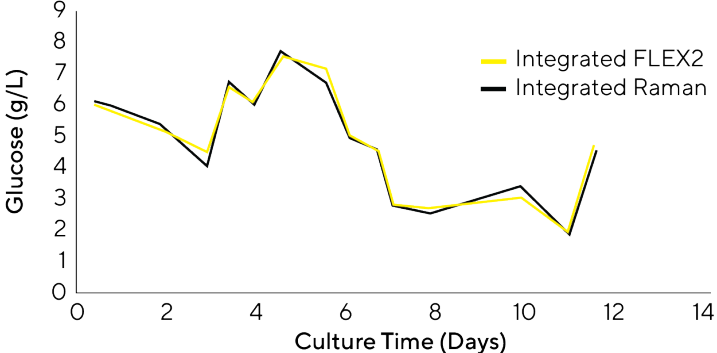
# Automated Feed Control in Ambr<sup>®</sup> Using Raman Spectroscopy



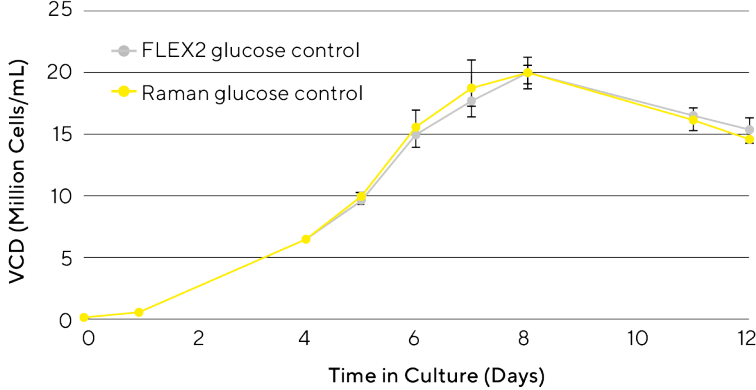
Glucose model built without spiking range: 2 – 9 g/L



Glucose model built with spiking range: 2 – 15 g/L



Raman predictions and NovaFlex 2 measurements give comparable results



Cell culture profiles are very similar for glucose control based on Raman or Nova FLEX2 assays

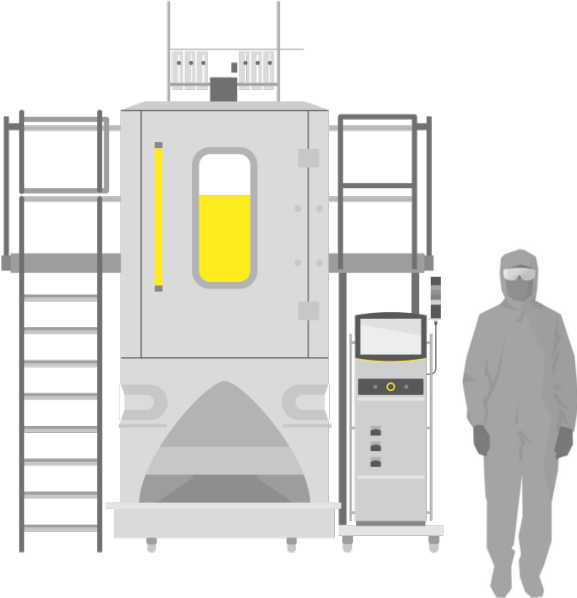
Data source: Ambr<sup>®</sup> 250 High Throughput, with Kaiser Raman spectrometer and probe; Marek Hoehse and Mike Sibley, Sartorius; Glucose set point: 5 g/L using bolus additions

# Raman Spectroscopy Use Cases

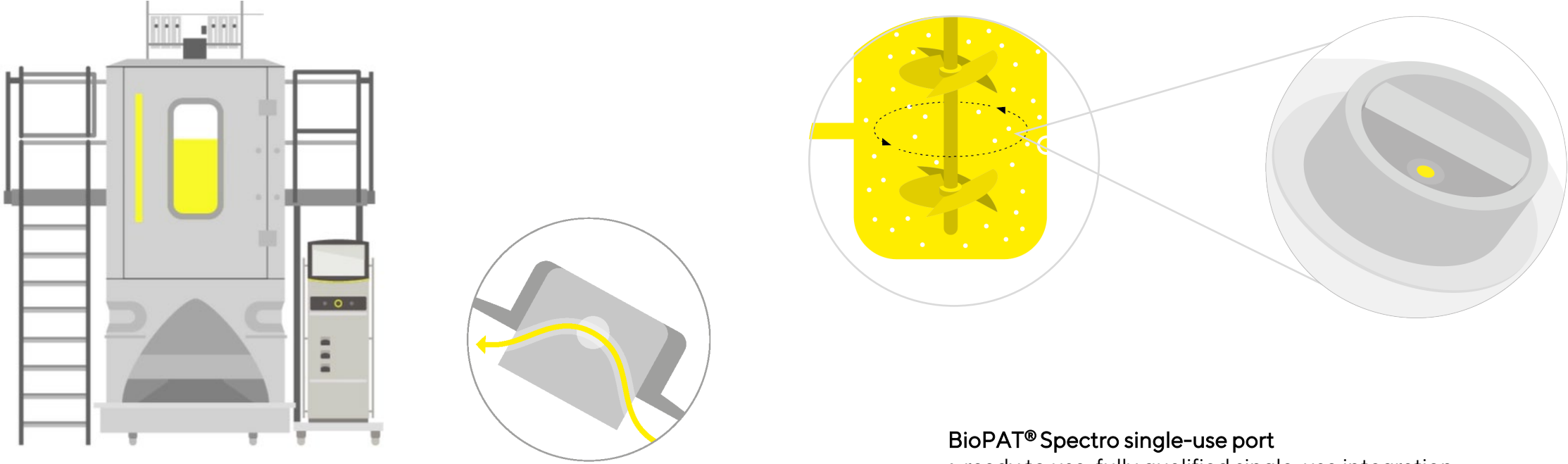
Model building

Process development

Single-use manufacturing



# BioPAT® Spectro Single-Use, Inline Integration in Flexsafe STR®



- BioPAT® Spectro single-use port**
- ready to use, fully qualified single-use integration
  - isolation of sample from light and bubbles
  - faster set-up time

# BioPAT® Spectro Single-Use, Inline Integration in Flexsafe STR®



BioPAT® Spectro single-use port in Flexsafe STR® bag



Connection of a Endress+Hauser Raman spectrometer probe



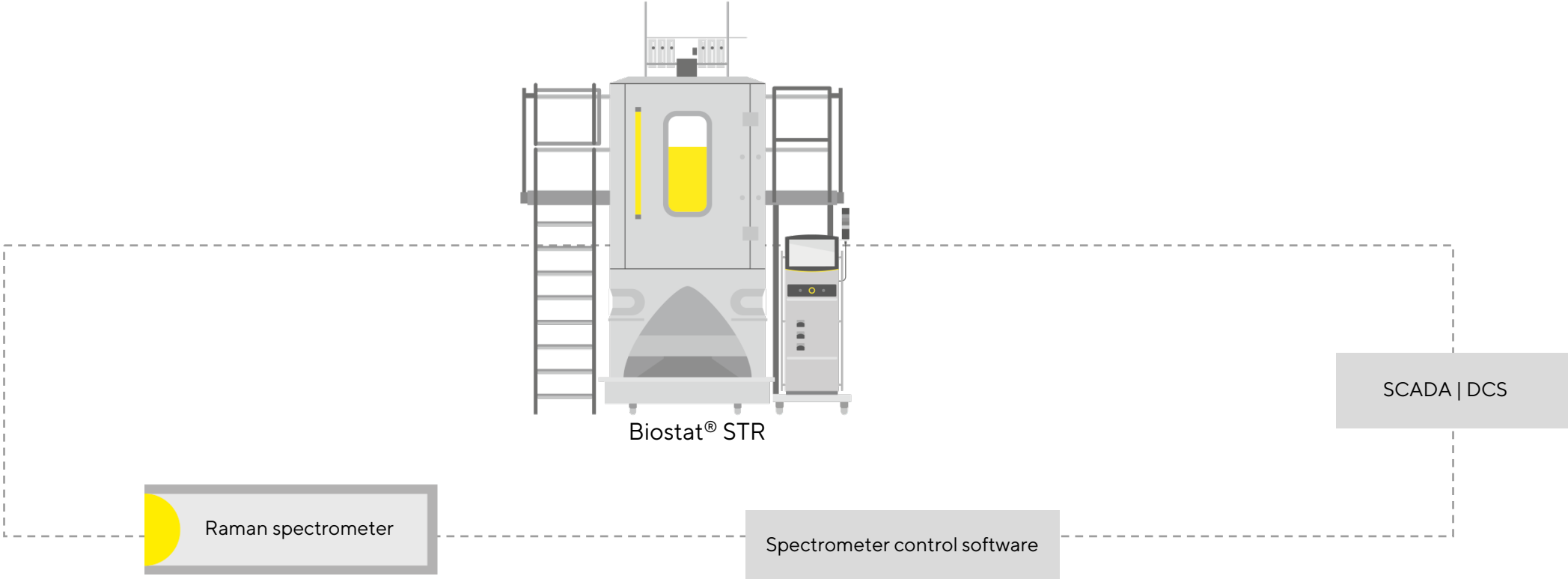
Connection of a Tornado Raman spectrometer probe)



Standardized optical design: same probe for Ambr® and Biostat STR®



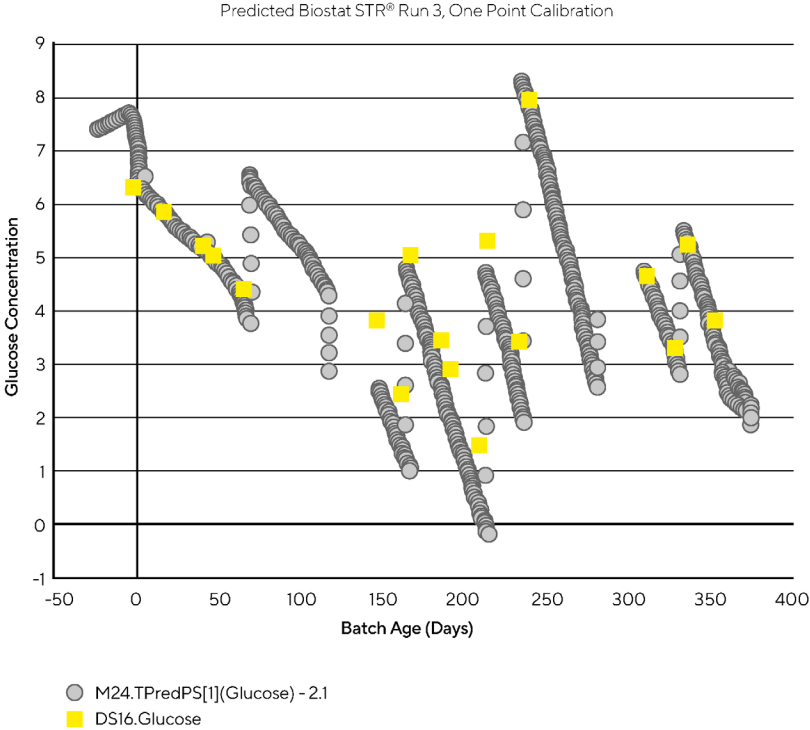
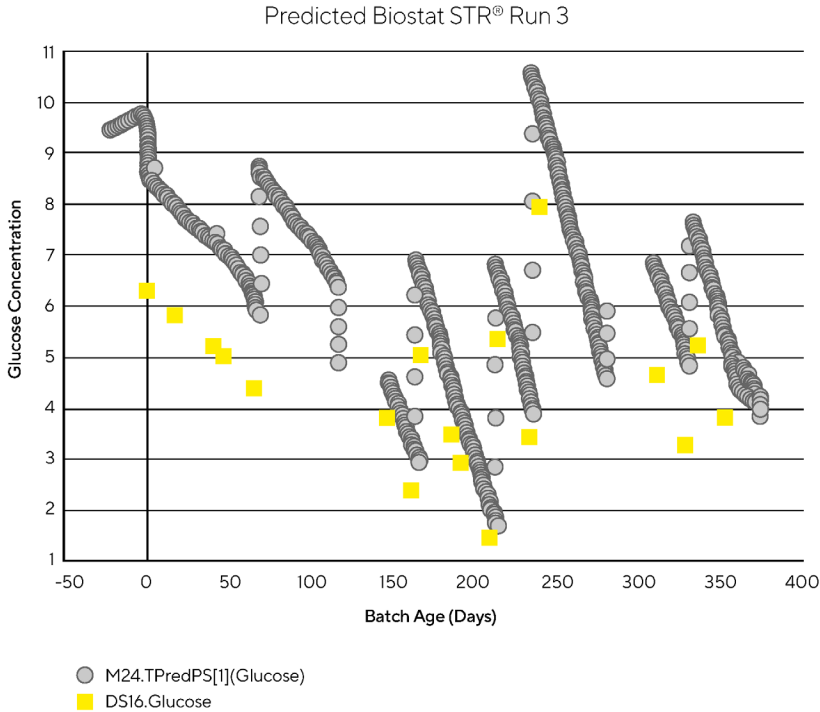
# BioPAT<sup>®</sup> Spectro integration in Biostat STR<sup>®</sup>



# Scalable Models Due to Platform Approach - Model Transfer to 50 L Single-Use Bioreactor

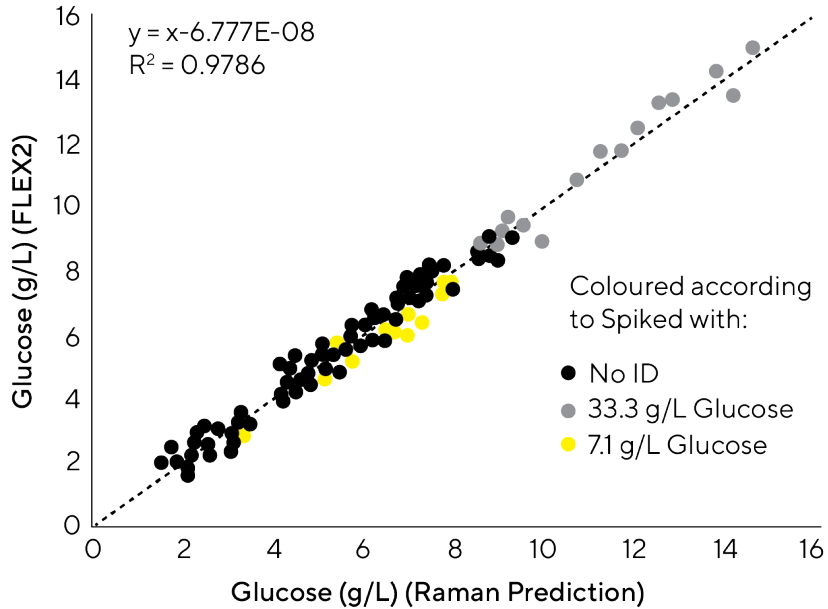
Model developed using Ambr<sup>®</sup> 15 runs + SUB batch 1  
Used to predict SUB batch 3

Potential for Model Transfer



Reproduced with kind permission from Ruth Rowland-Jones, GSK Stevenage. Performed with a prototype spectroscopy integration and a Tornado Raman spectrometer.

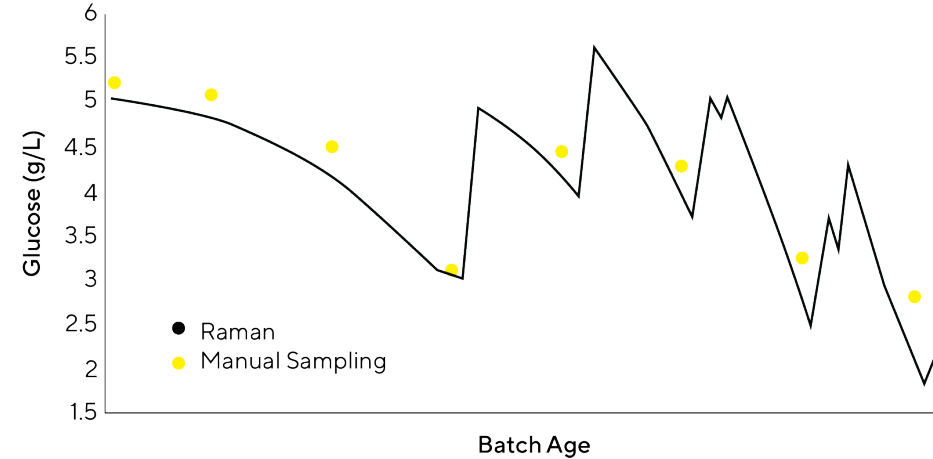
# The Platform Approach Enables Model Transfer Across Scales



## Model building in Ambr® 250

- Glucose model built using two Ambr® runs of eight vessels each
- Spiking two glucose stock solutions
- RMSE<sub>cv</sub> = 0.34 g/L

Data source: Ambr® 250 High Throughput & Flexsafe STR® 200, with Kaiser Raman spectrometer and prototype probe; Marek Hoehse and Mike Sibley, Sartorius



## Direct model transfer from Ambr® to Flexsafe STR® 200L

- STR and Ambr® runs happened at different sites
- Different seed train, media lot and reference method were used
- The raw Ambr® model was used without any additional data from an STR
- Offset correction was performed based on first reference measurement

# Enable Rapid and Robust Scale-Up With Scalable Bioreactor Hardware

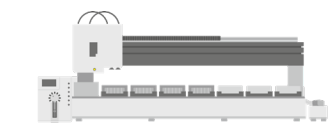
- From Ambr® 250 to Biostat STR® 2000 Sartorius vessels maintain geometric similarity
- Ensures that mixing and oxygen transfer rates remain consistent over scaling
- Simplifies technology transfer between cultivation scales



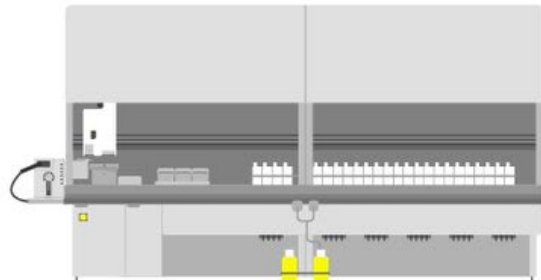
Ratio	Ambr® 250	Biostat STR® 50	Biostat STR® 200	Biostat STR® 500	Biostat STR® 1000	Biostat STR® 2000
Vessel Height/Diameter	2	1.8	1.8	1.8	1.8	1.8
Impeller Diameter/Vessel Diameter	0.42	0.38	0.38	0.38	0.38	0.38

# BioPAT<sup>®</sup> Spectro: Scalable Raman Spectroscopy in Single-Use

- Robust model building in Ambr<sup>®</sup> is faster, easier, and more cost effective than possible before
- Quality by design is enabled by introducing Raman spectroscopy to high-throughput process development
- Efficient transfer to single-use manufacturing due to the scalable platform approach



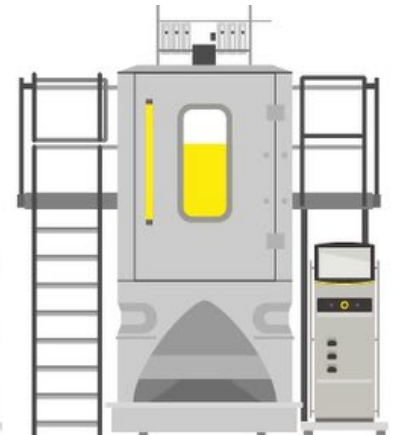
Ambr<sup>®</sup> 15 Cell Culture



Ambr<sup>®</sup> 250 High Throughput



Biostat<sup>®</sup> STR



# How to Get BioPAT<sup>®</sup> Spectro?

- All Ambr<sup>®</sup> 250 High Throughput and Ambr<sup>®</sup> 15 Cell Culture instruments can be retrofitted
- Quotes can be generated right now, orders will ship starting from 08/2020
- Flexsafe STR<sup>®</sup> bags with the BioPAT<sup>®</sup> Spectro single-use ports are available from 12/2020 onwards
- Please contact your local PAT expert for inquiries

<b>Name</b>	<b>Region</b>	<b>Contact</b>
Michael Isaacs	North America	Michael.Isaacs@Sartorius.com
Sabah Shaker	EMEA	Sabah.Shaker@Sartorius.com
Jiahui Yu	Asia	Jiahui.Yu@sartorius.com

# Thank you.

**Veronica Brivio**  
Product Manager PAT  
Mobile ++44.1763.2274.88  
[Veronica.Brivio@Sartorius.com](mailto:Veronica.Brivio@Sartorius.com)

**Dan Kopec**  
PAT Expert North America  
Mobile +1.480.251.9049  
[Dan.Kopec@Sartorius.com](mailto:Dan.Kopec@Sartorius.com)

The Sartorius logo is displayed in a bold, black, sans-serif font on a bright yellow background. The letters are closely spaced and have a clean, modern appearance.