

# SARTORIUS

## Simplifying Progress

### OSIsoft and Sartorius Partner to Help CDMOs Leverage their Data

November 24th, 2020

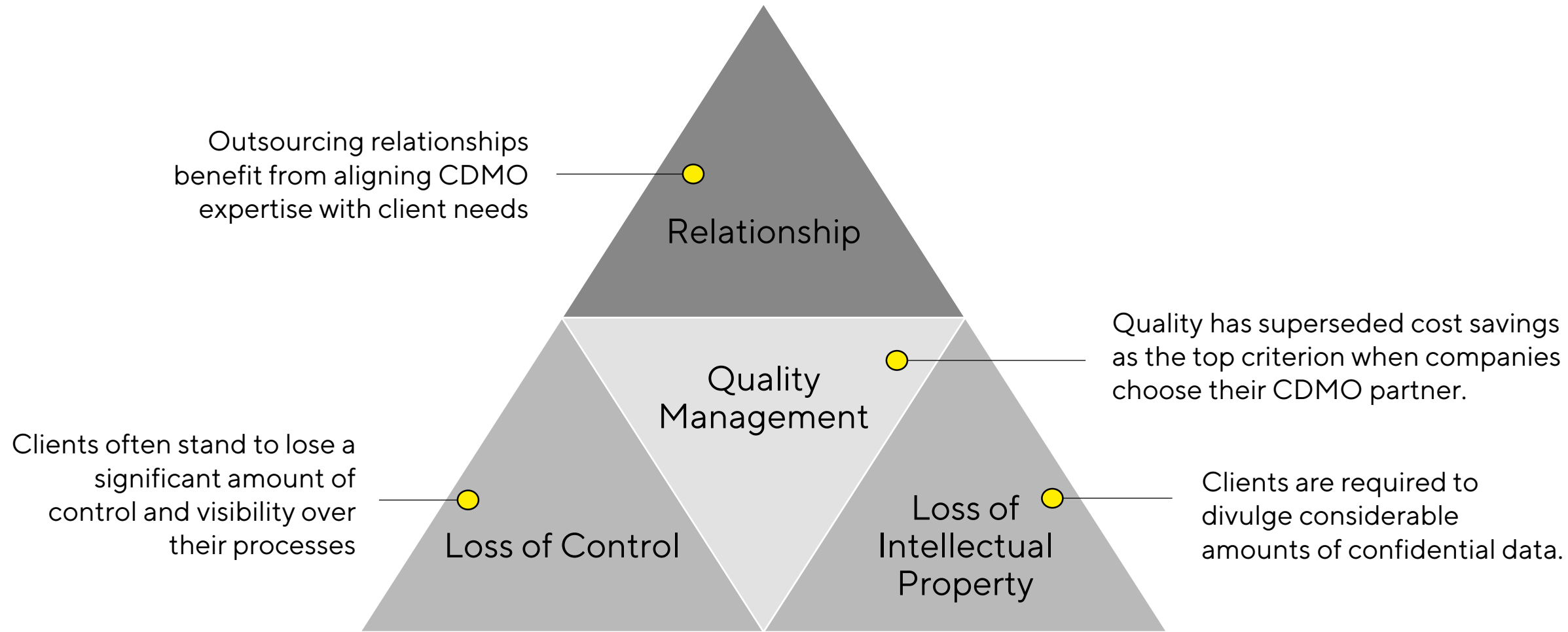


# Agenda

- Introduction: Data integrity: The New CDMO Challenge
- Presentation from OSIsoft
  - Who is OSIsoft?
  - OSIsoft in Life Science Industry
  - How to handle Data Integrity in a Global Ecosystem
  - OSIsoft PI System Implementations at CDMOs
- Faster and More Reliable Operations: A FUJIFILM Diosynth Biotechnologies' Case Study
- Closing: Q&A



# If you Remember from our Last Webinar...



# New CDMO Challenge: Data Access and Sharing

Disparate internal and external IT and data management systems

Data stored in non-electronic systems (e.g. paper records)

Increased complexity in global supply chains

Compliance in a new era of regulatory scrutiny (FDA & SEC)

Process complexity of new therapies



Top Data Related Challenges in the CDMO Industry



# So What?



These CDMOs struggle to provide secure, compliant data sharing with their partners



# Lack of Data Transparency Can Result In....

- **Delay in Business Critical Processes** – tech transfer, IP transfer, regulatory approval,...
- **Operational Inefficiencies** – longer cycle times, lower yields, minimized throughput...
- **Cost to Reputation** – failed audits, extended time-to-market, supply chain shortages...





It's your data - take control

WHO WE ARE

# Pioneers in operational data management for essential industries



**40**

years

**\$5B**

valuation

**21k**

sites

**140+**

countries

**2B**

data streams

**24\***

of our first 25 customers  
are still with us



# Leading the market in critical operations



## Oil & Gas

**85%**

of the top oil and gas companies



## Power & Utilities

**1000+**

utilities worldwide



## Mining & Metals

**25**

of the top 25 mining and metal companies



## Pharma & Life Sciences

**25**

of the top 25 pharmaceutical companies



## Chemicals

**9**

of the top 10 chemical companies



## Manufacturing

**6**

of the World Economic Forum's Factories of the Future

# OSIsoft: Proud to support the Pharmaceutical Industry

Frame = EA customer



# OSIsoft: Proud to support the Pharmaceutical Industry



ALFASIGMA

Perrigo

Alkermes

GCBT B|BRAUN

Ortho Clinical Diagnostics  
a Johnson & Johnson company

stryker

FERRING  
PHARMACEUTICALS

InnoTech  
ALBERTA

WuXi Biologics  
Global Solution Provider  
A WuXi AppTec Group Company

AUDENTES  
THERAPEUTICS

Ampio  
PHARMACEUTICALS

VIR

Intarcia  
Therapeutics, Inc.

BIO-RAD

PAREXEL

INTERQUIM

Recipharm  
good for business

HUMACYTE

FUJIFILM  
Diosynth  
biotechnologies

MILLENNIUM PHARMACEUTICALS, INC.

brammer  
Manufacturing Personalized™

VERTEX

Paragon  
BIOSCIENCES

MACROGENICS

lek  
a Sandoz company

FMC ZENTIVA

GEDEON RICHTER

POWERED BY SCIENCE  
GENENCOR

天士力 TASYL

Alcon

SK biotek

KBI  
BIOPHARMA

West

BioVectra

Seqirus  
A CSL Company

DePuy Synthes  
COMPANIES OF Johnson & Johnson

bluebirdbio

# Digital Fundament

Plug & Produce



Knowledge driven



User Driven

Agile  
Transparent  
Value driven



Information

100  
010

Data centric

- Accesible
- Scalable

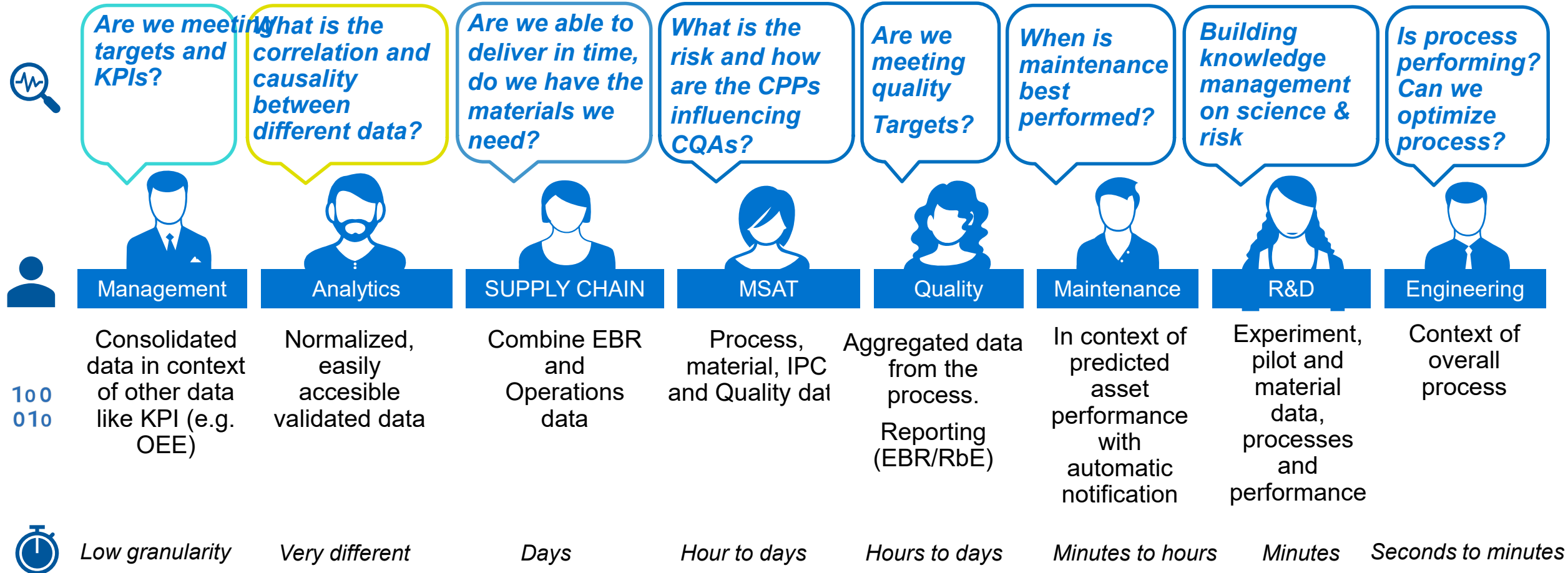


Metadata

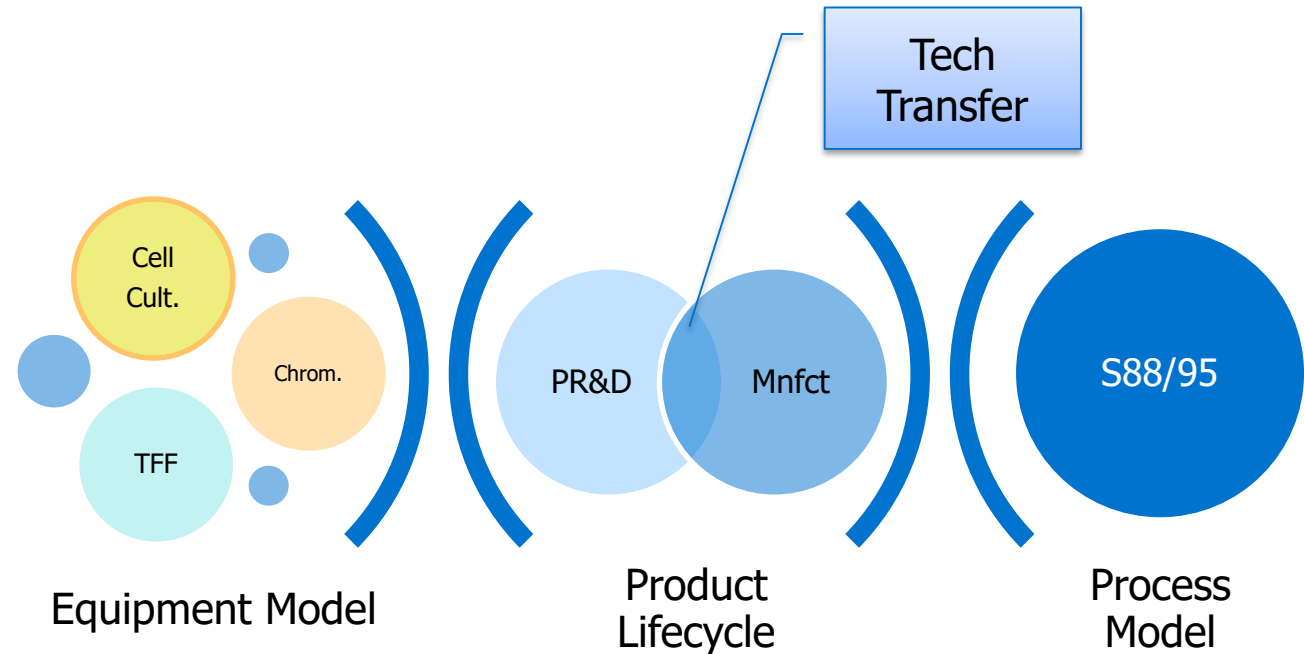
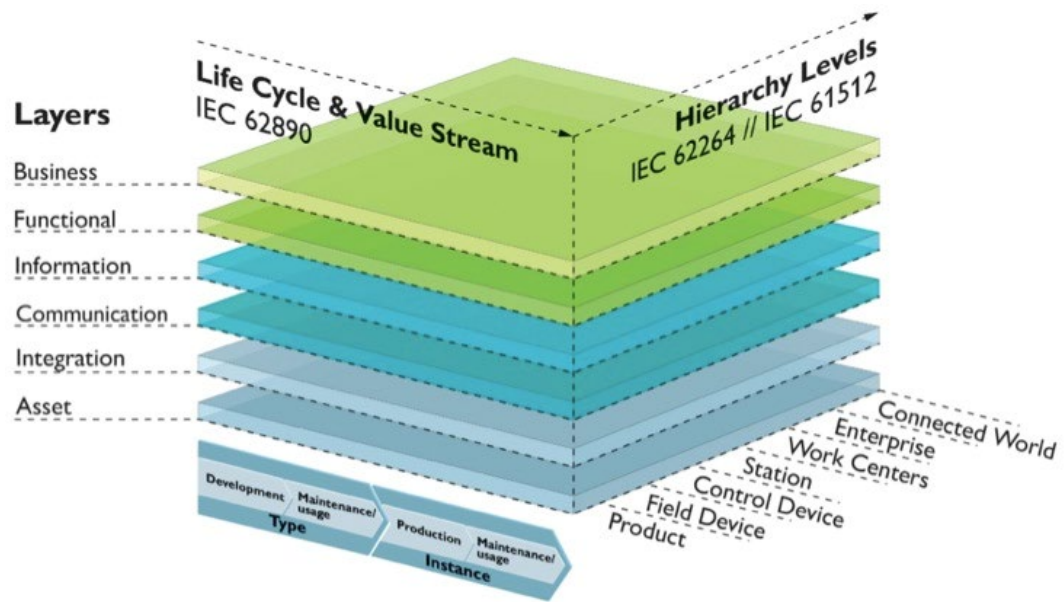


Data

# User Driven, who are the drivers?



# Three key dimensions



# Equipment & Process models builds the Smart

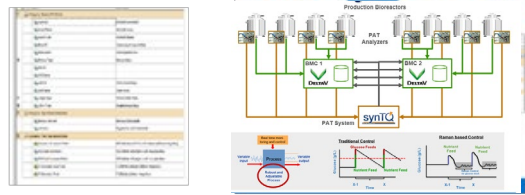
## Pharma



Fermentation Template



Chrom. Template



Smart Application Template

- Metadata
- Sensors
- Time Series
- Equipment data
- Performance curves
- KPI / OEE
- SOP
- Alarms
- Events
- Notifications
- Predictions
- Biomass
- Product
- Metabolites
- Endpoints
- Calibration status
- Etc.

**E.g. Bioreactor Digital Twin**



Bio Pharma Plant #1



Pharma Plant #2



CDMO #1

- Ozona
- Acid gas Removal
- Dehydration
- Fractionation Train
- Gathering Systems
- Mercury Removal
- Meter stations
- Nitrogen Reinjection
- Sulfur Unit
- Sweetening Unit
- Tail gas Treating

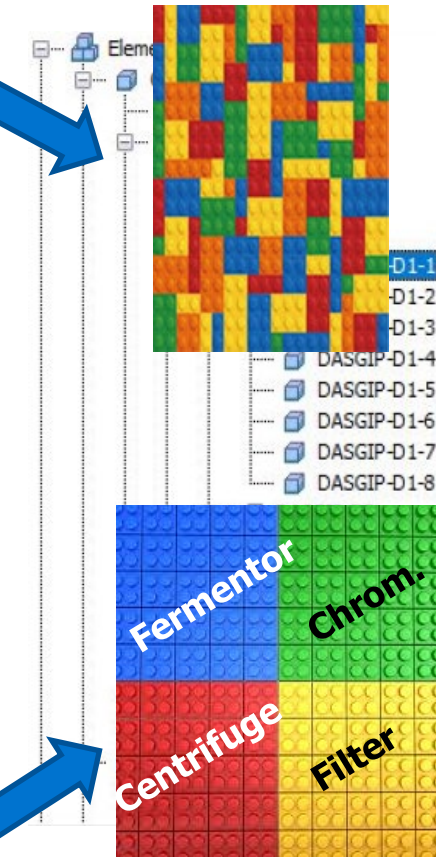
Digital Plant #1

- Ozona
- Acid gas Removal
- Dehydration
- Fractionation Train
- Gathering Systems
- Mercury Removal
- Meter stations
- Nitrogen Reinjection
- Sulfur Unit
- Sweetening Unit
- Tail gas Treating

Digital Plant #2

- Ozona
- Acid gas Removal
- Dehydration
- Fractionation Train
- Gathering Systems
- Mercury Removal
- Meter stations
- Nitrogen Reinjection
- Sulfur Unit
- Sweetening Unit
- Tail gas Treating

CDMO #1



Digital Enterprise

# Pharma, a truly global ecosystem

-  Key Data Centre
-  Manufacturing Plant
-  CMO/CDMO/CMA/CRO

**The Numbers Game**

- > 200 Individual Product Families
- > 2000 SKUs
- > 18000 Raw Material SKUs
- 29 Internal Sites
- 130 External Sites
- 0 Products Made all Internally

	External Mfg	Procurement	Total
# of Partners	108	480	588
# of Sites	130	614	744
# of SKUs	2000	18,000	20,000

janssen



# Global Process Maps



## Process Visualisation



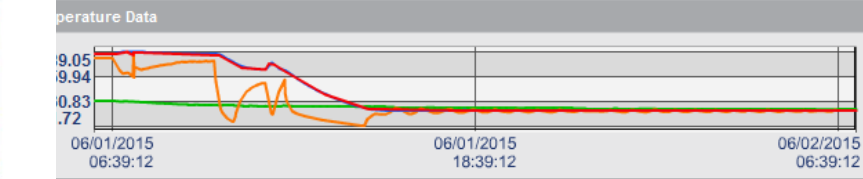
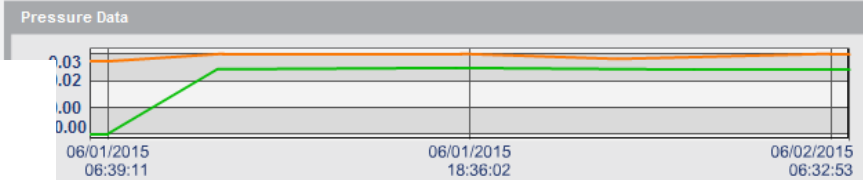
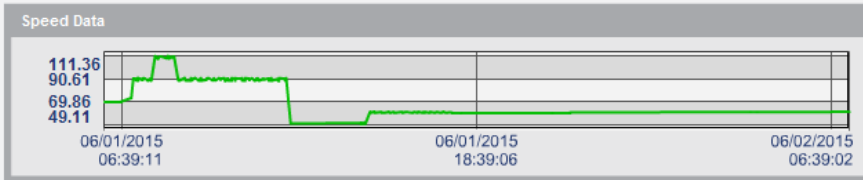
Unit Selection

Cork Chemical: Reactor - R221

Schedule Attainment

Detailed Analysis

All Sites  
 All Sites  
 Biogen Idec  
 Beerse  
 Cork Bio  
 Cork Chemical  
 Plant 3  
 Mod1



FOR INFORMATION ONLY, NOT FOR GMP USE.

Speed (RPM)	60.36	Low Pressure (psi)	0.03
High Pressure (psi)	0.03	Jacket Temp (°C)	19.85
Vapour Temp (null)	20.97	Middle Temp (°C)	19.79
		Bottom Temp (°C)	20.07

Global Interconnected Standard Visualization

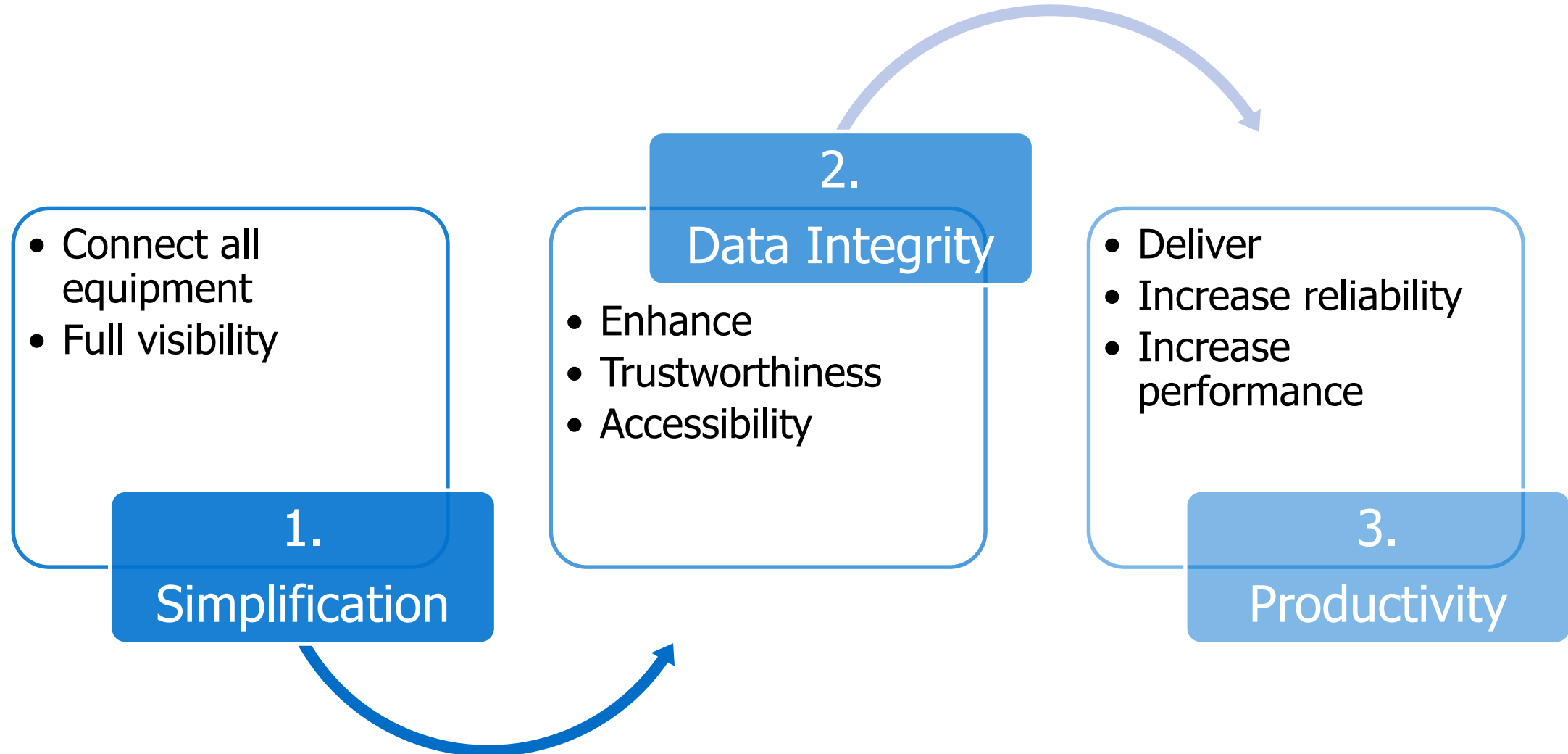
One  
PI SYSTEM

Source: P McKenzie Janssen

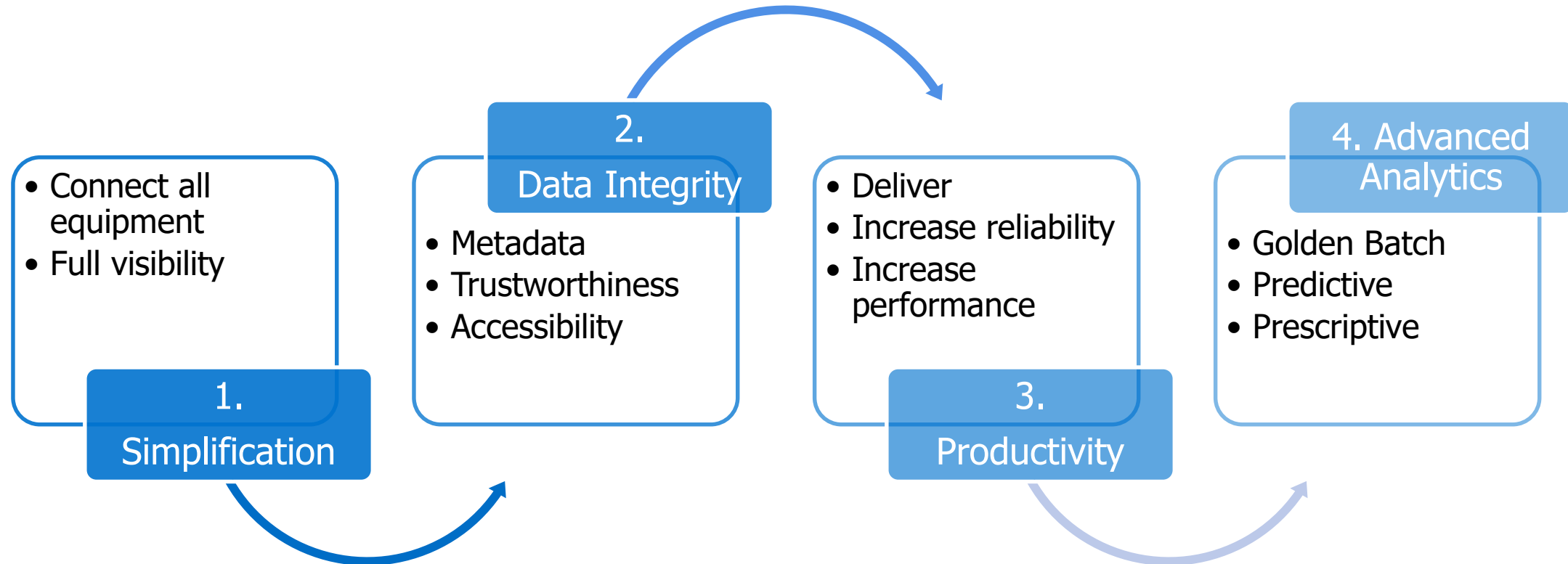
# How do we reach there?

Remember it is about the business strategy

# From standalone to integrated operation

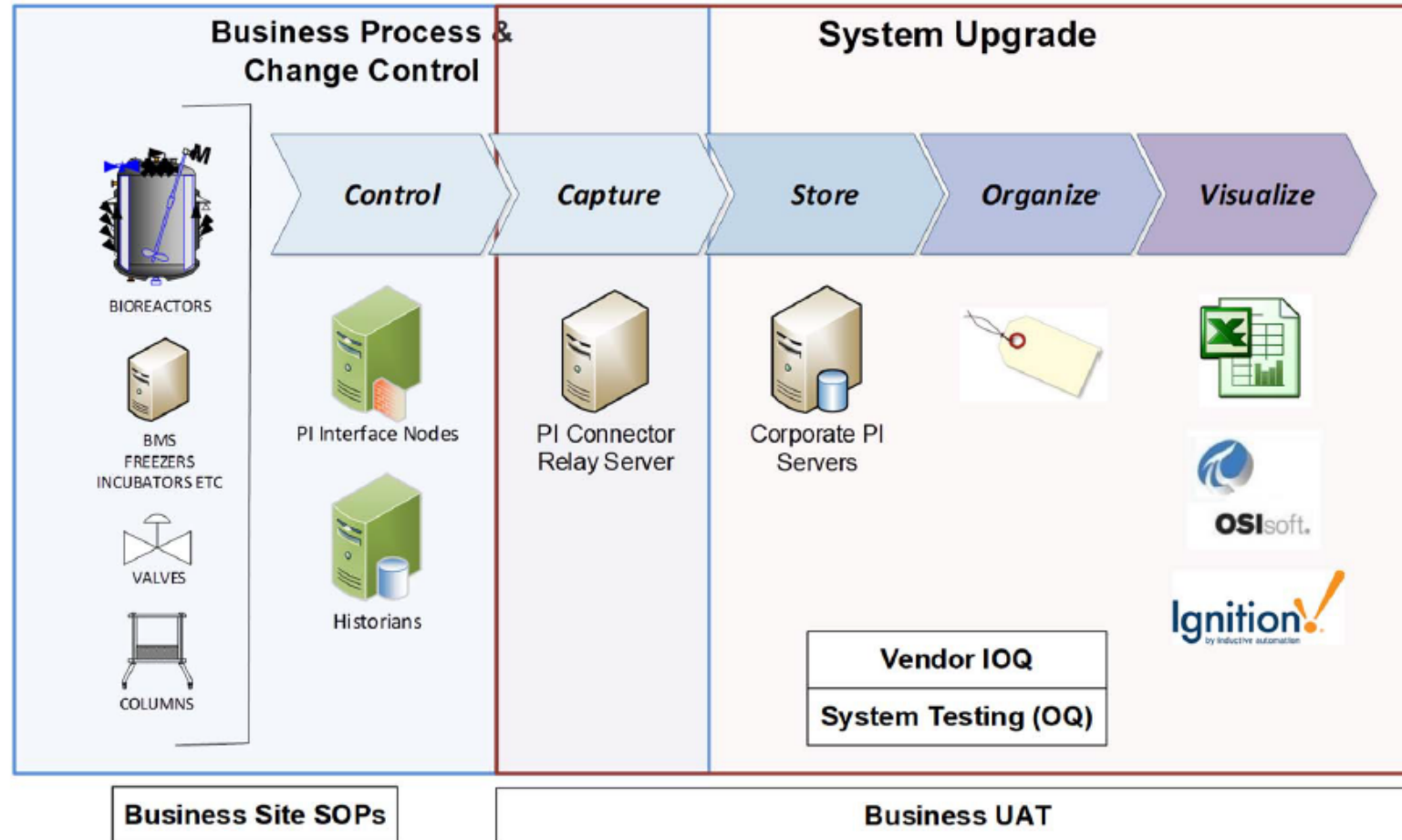


# From standalone to integrated operation



# 1. Simplification

# PI Enterprise Validation Approach

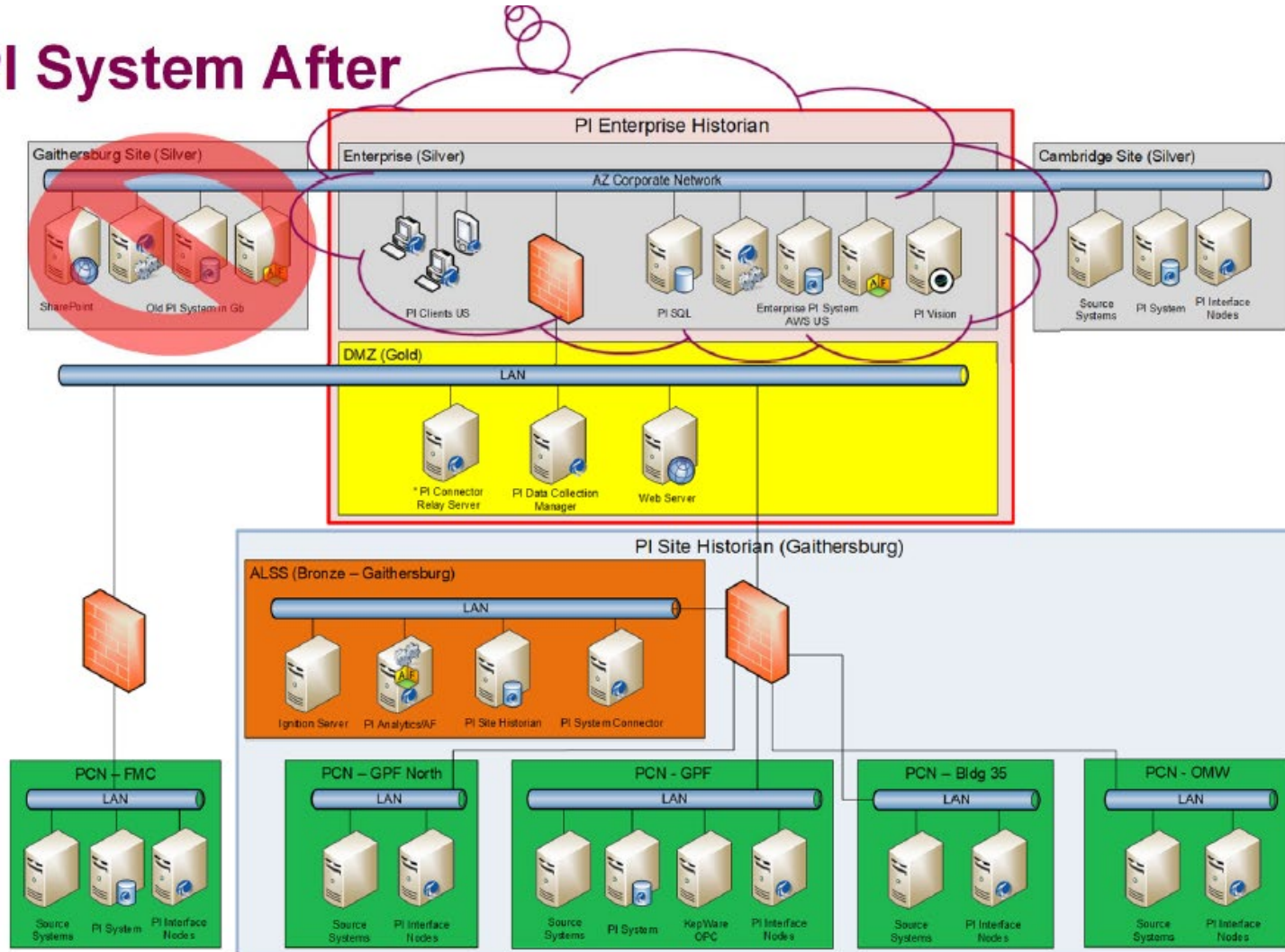


15

❖ Focus on the Enterprise Historian NOT the Site

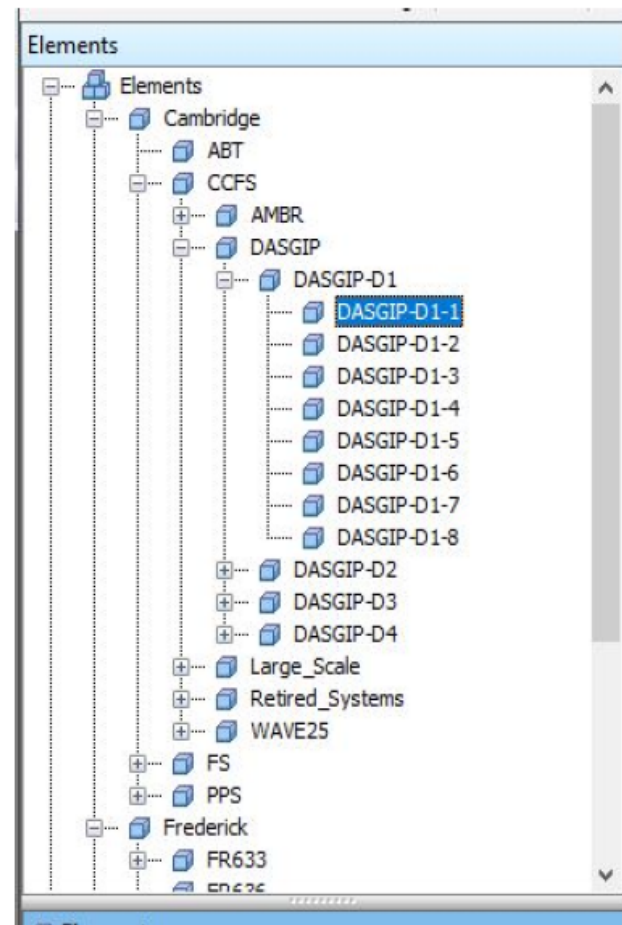


# OSI PI System After



## Customer Value – Tags & Assets

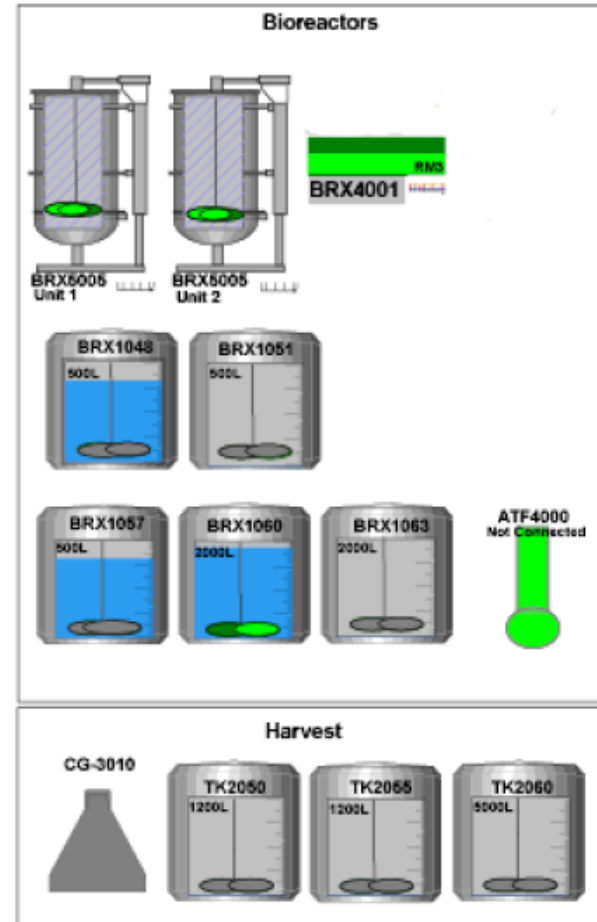
- ❖ Deployment of new assets with templates
- ❖ Standardized navigation for assets allowing users to quickly find instruments and data of interest
- ❖ Ability to integrate process data with other asset context
- ❖ Alignment with AZ standards for tag naming and architecture





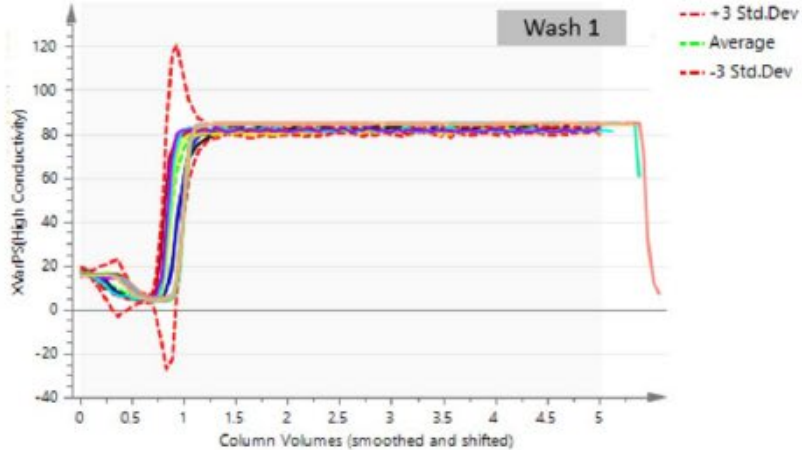
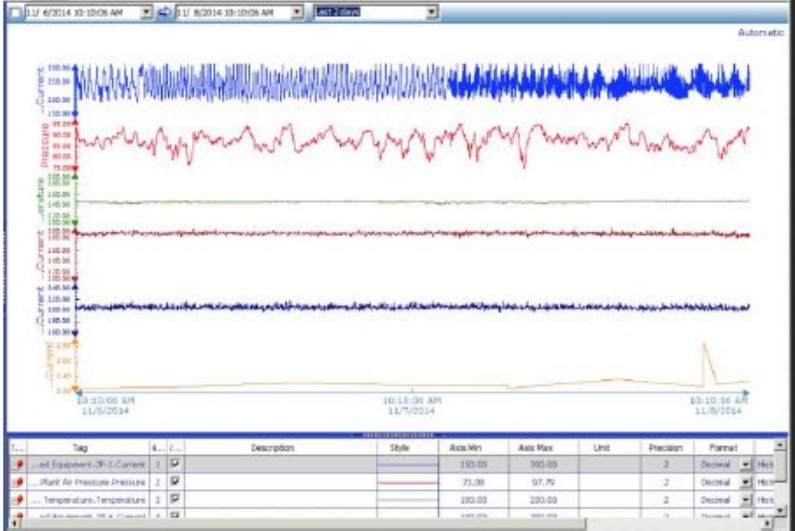
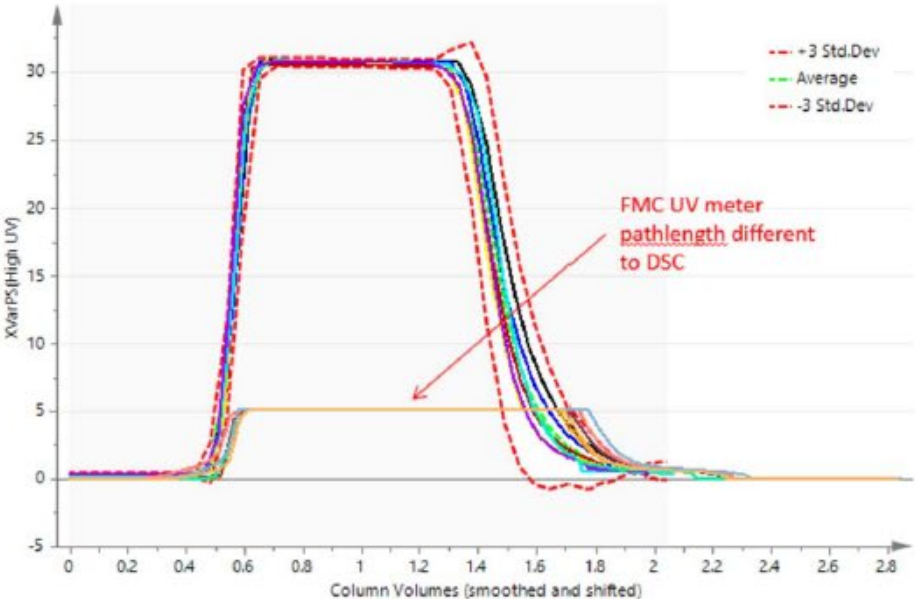
# Customer Value – User Interface

- ❖ An enhanced, web-based, user experience
- ❖ Increased application stability
- ❖ Ad hoc report creation without needing local PC installs, reducing overhead for our support teams and our users
- ❖ Improved reporting for GMP equipment
- ❖ No interruptions to critical processes in our R&D labs and for pilot plant operations



# Customer Value – Future

- ❖ AWS cloud supports better collaboration across sites
- ❖ Capability to determine “Golden Batch”
- ❖ Ability to use mobile devices



# 2. Data Integrity

# Validation of a global data mart

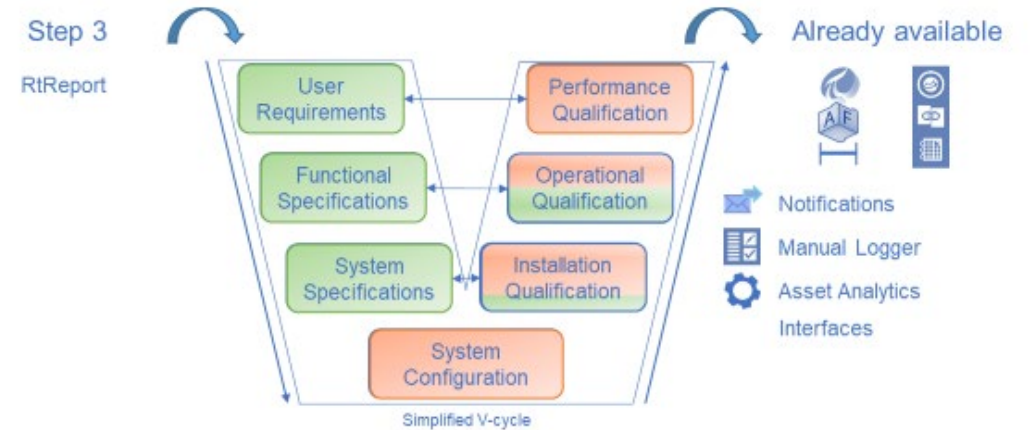
## GlaxoSmithKline Vaccines

Creating a global PI Solutions Store to Ease Site Deployment and Increase Compliance

CHALLENGE	SOLUTION	RESULTS
<p>Align and accelerate OSIsoft PI deployment in parallel on several manufacturing sites</p> <ul style="list-style-type: none"> <li>Several sites implementation to manage in parallel</li> <li>Minimize the local efforts for implementation &amp; qualification</li> </ul>	<p>Global qualification approach of OSIsoft tools with pre-qualified and documented modules</p> <ul style="list-style-type: none"> <li>Strong partnership with OSIsoft (EA) and PAS integrator</li> <li>Involvement of local sites in the budget estimate and scope definition</li> </ul>	<p>60% reduction of local qualification effort</p> <ul style="list-style-type: none"> <li>Similar approach for other systems: ROI after 5 instances implementation</li> <li>Expected to quickly increase the solutions available thanks to the collaborative approach</li> </ul>



## The infrastructure qualification approach



### Solutions

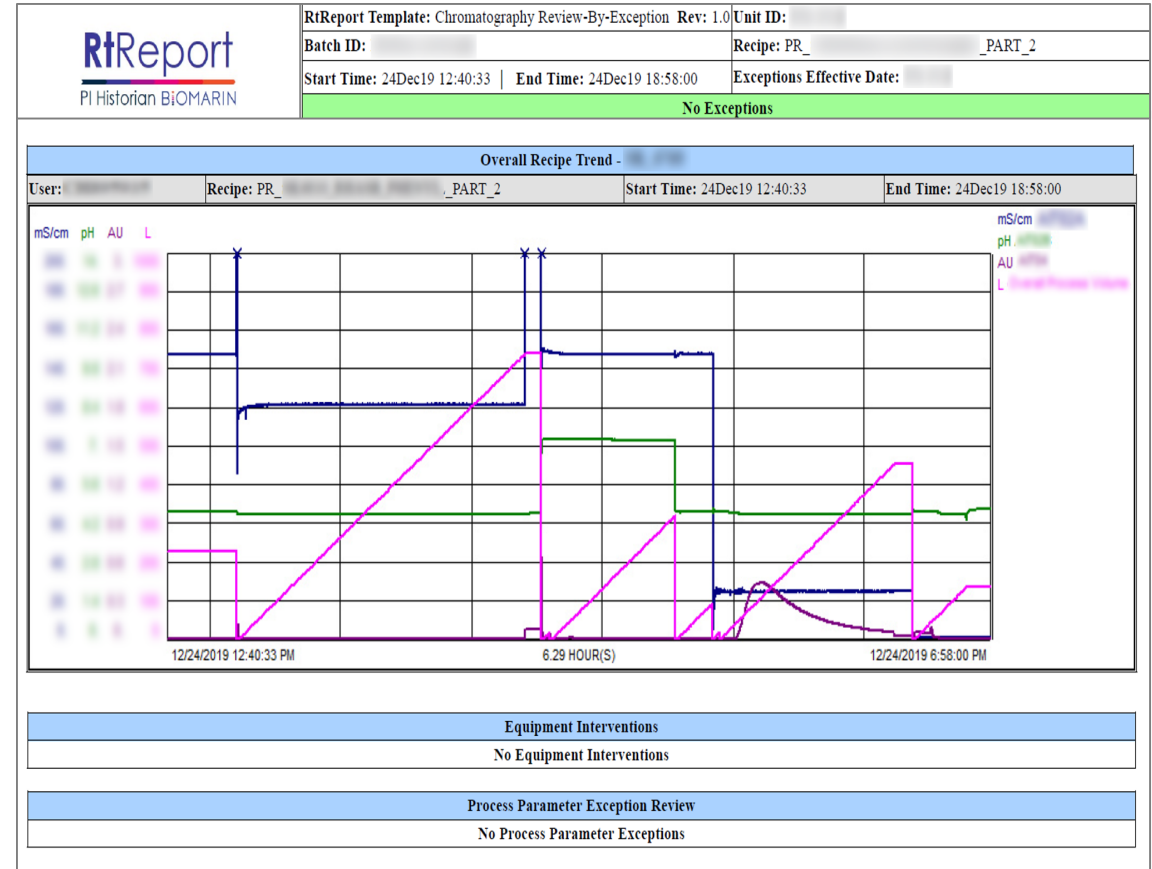
<b>Facilities</b>	<b>Maintenance 4.0</b> CBM & Predictive	<b>Alarms Analytics</b> Equipment & PCS	<b>EMS Analytics</b> Room analytics / HVAC Qualification
	<b>Deviation Investigation</b> Pi-DataLink, Pi-Vision & RT-Reports	<b>Energy Analytics</b>	
<b>Equipments (Std objects)</b>	<b>Autoclaves</b>	<b>Fermentors</b>	<b>Filling Line</b>
	<b>Reporting</b> (GMP or not)	<b>Digital Dashboard</b> Realtime data display	<b>Process Analytics</b>
	<b>Machine Learning</b> OEE, Analytics ...		

# Data Integrity

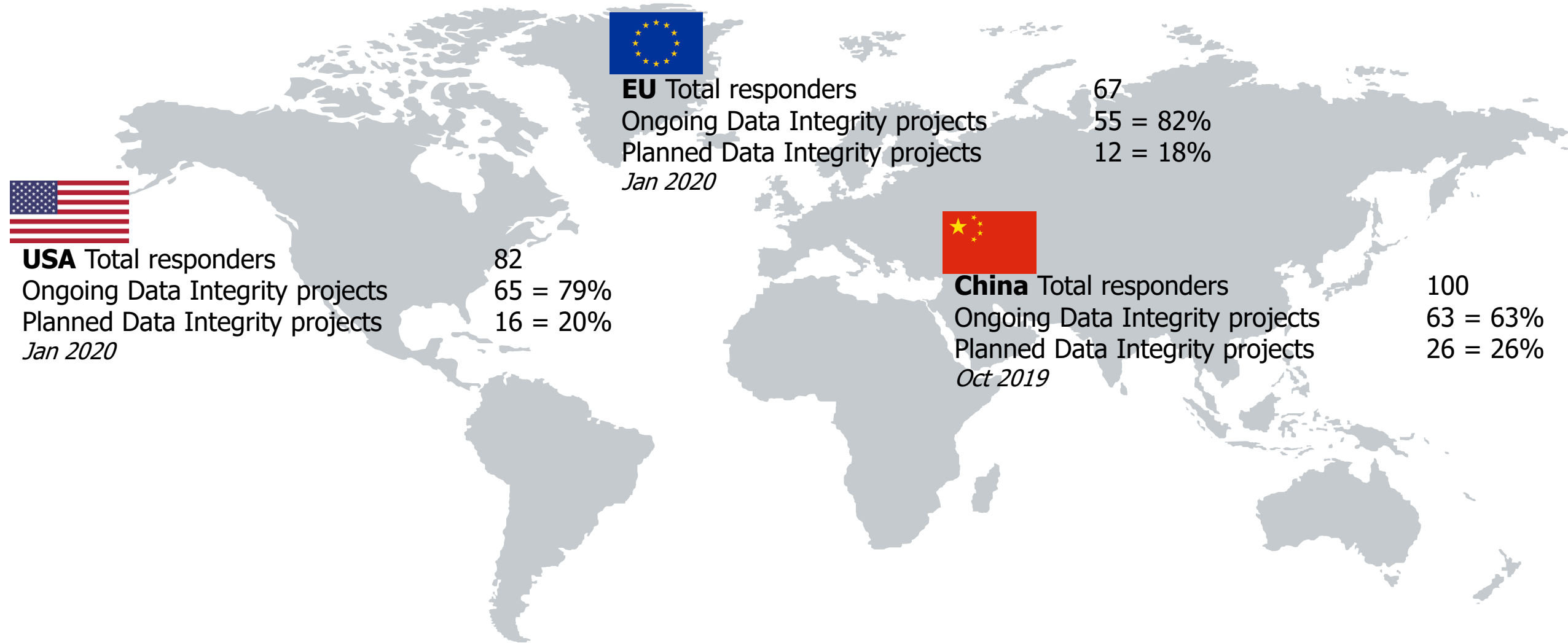
## Before - Paper



## After - RbE



# Data Integrity – a global snapshot



# Data Integrity and Contract Organizations (CMO/CDMO)

**Carmelo Rosa, Director of FDA OMPQ's**

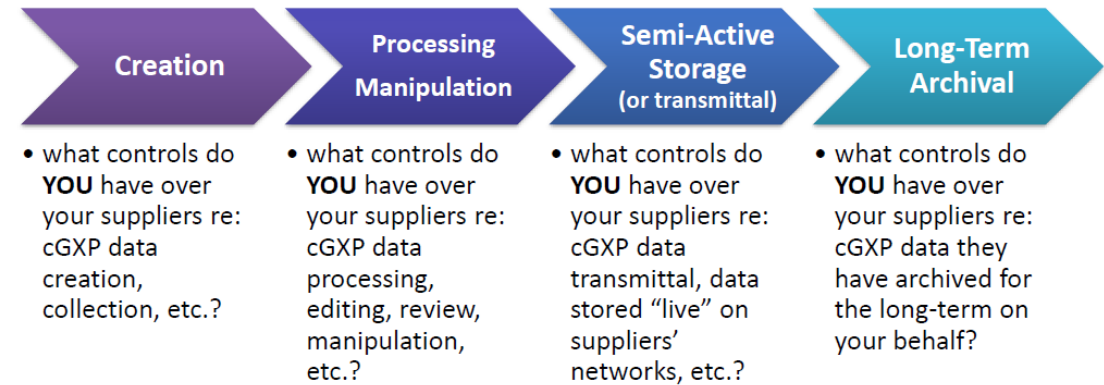
*"Data integrity issues have always existed!"*



Drug makers should not look to contract manufacturers to reduce their responsibility for data accuracy and reliability, Some biopharma companies regard contract testing and production operations as one way to alleviate their involvement in inspections and dealings with regulatory authorities.

**Data Integrity issues are a Global problem**

## Data Integrity Lifecycle



© 2016 Cerulean Associates LLC

www.ceruleanllc.com

8

Source: **John Avellanet – CMO Conference** [www.ceruleanllc.com](http://www.ceruleanllc.com)

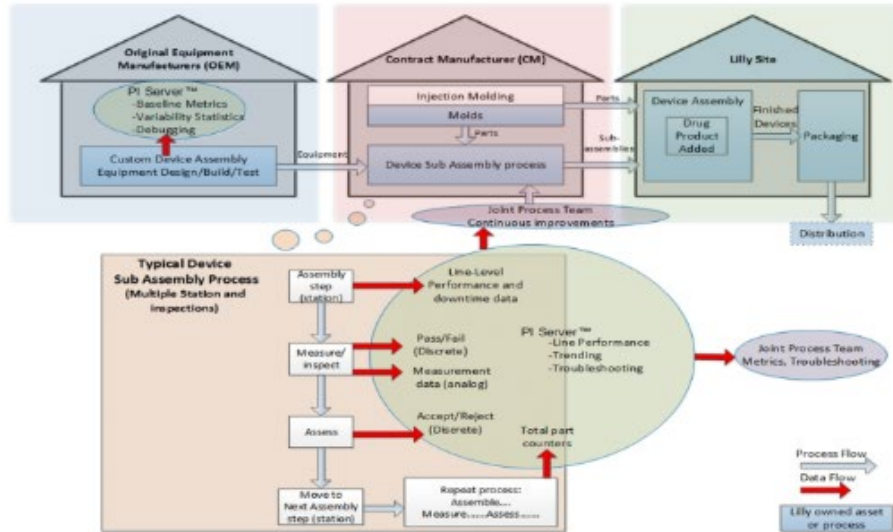
*"Every gram we produce is valuable, however the data behind every gram is more valuable"* SVP Lonza

# 3. Productivity



# Supply chain management and CMO

## OEM to CMO to Lilly – Supply Chain



Real-time visualization of the process data in a format that's easy to understand from the CMO.



# 总结和概览

**WuXi Biologics**  
Global Solution Provider

无锡药明生物技术股份有限公司

PI系统作为智能应用平台用于监控、分析和优化工厂的设备 and 工艺，提高批次工艺的稳定性，并将逐步应用于全球生产基地。

As an intelligent application platform, the PI system is used to monitor, analyze and optimize plant equipment and processes, improve the stability of batch processes, and will gradually be applied to global production bases.

刘松 自控与信息管理部 执行主任 - Liu Song Executive Director of Automation and Information Management Department



## 业务挑战

Continuous process data cannot provide batch analysis

The traditional preventive equipment maintenance program cannot reflect the actual operating status of the equipment, and it is easy to cause production losses when the equipment fails

There is no reference case for the construction of process system equipment and equipment model

## 解决方案

Build batch management of cell culture & purification and liquid preparation system

Use PI's AF and EF to generate event and operation records of equipment in the dosing system

Real-time monitoring of the operating conditions of the equipment in the operating cycle and early warning (including more than **6000** valves, **3000** gaskets, more than **100** pumps and agitators)

## 客户成效

Process batch management can better provide process monitoring and deviation warning

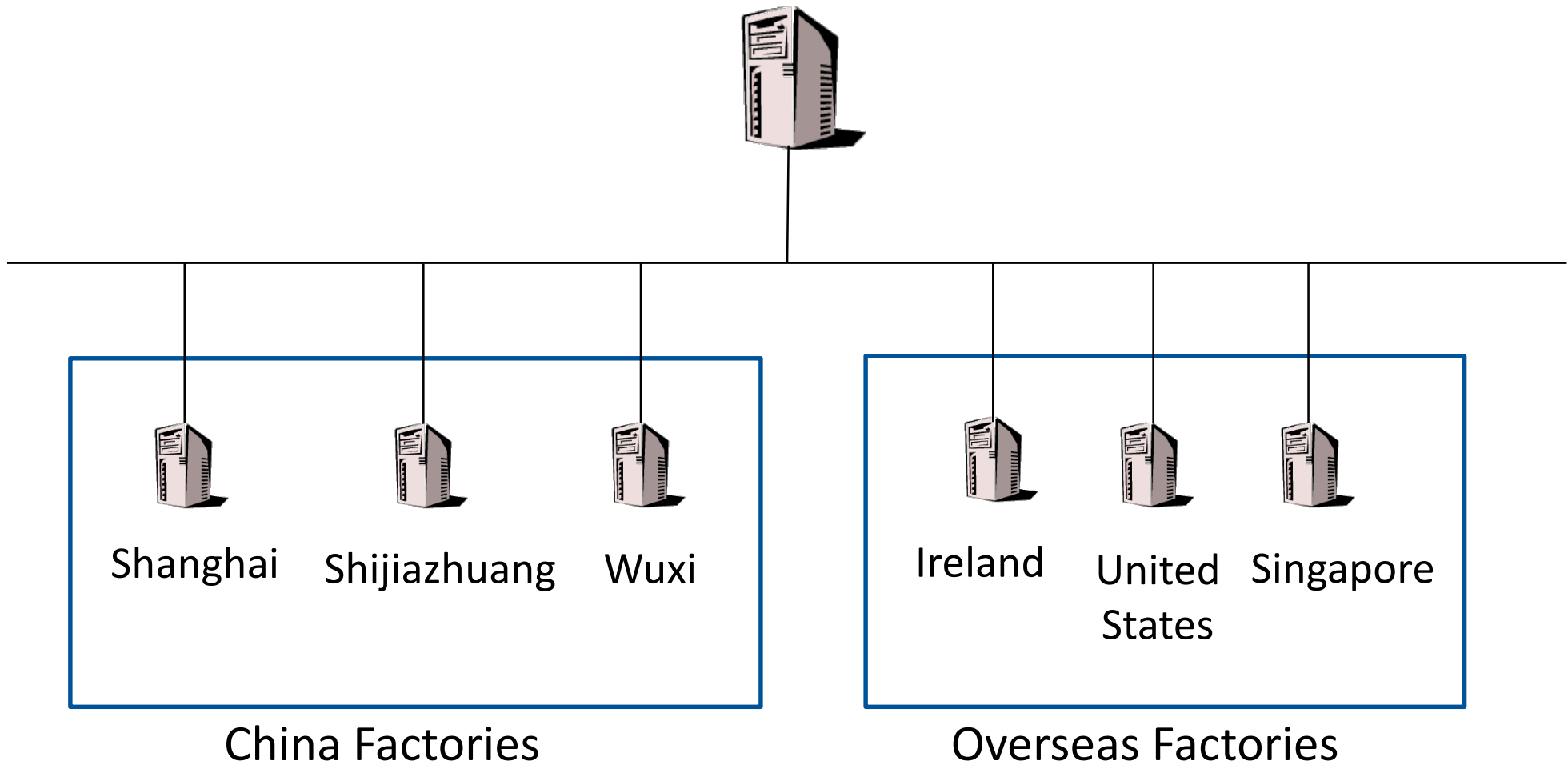
Monitor the health factor of equipment in real time and guide equipment maintenance based on early warning to reduce the cost of equipment maintenance

Transform preventive maintenance based on equipment usage to predictive maintenance based on equipment status



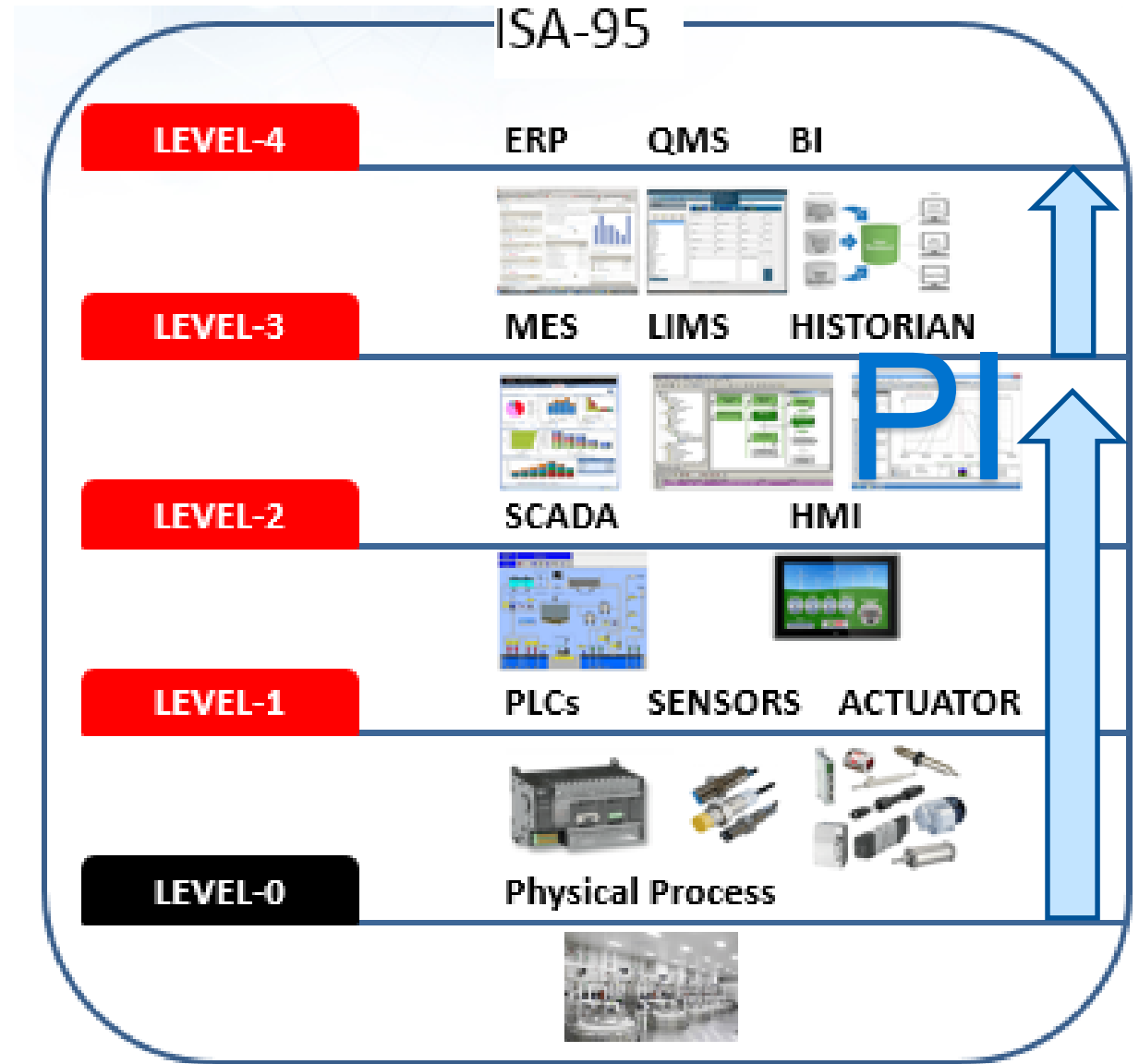
# PI System Global Architecture Diagram

Worldwide HQ ( Wuxi, China)



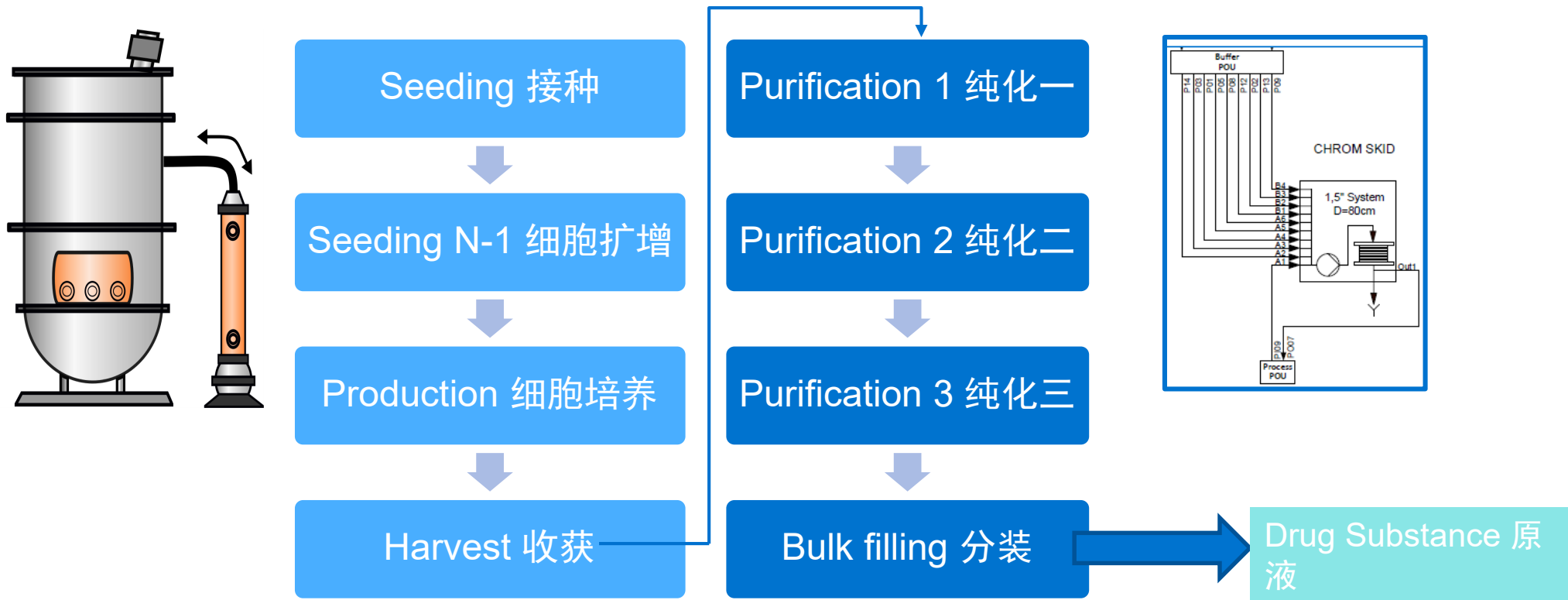
# PI 系统策略

- PI system is more than a historian, PI system is more than a reporting tool.
- A production data platform, provide data access, digital management and analysis
- GMP or Non-GMP ?

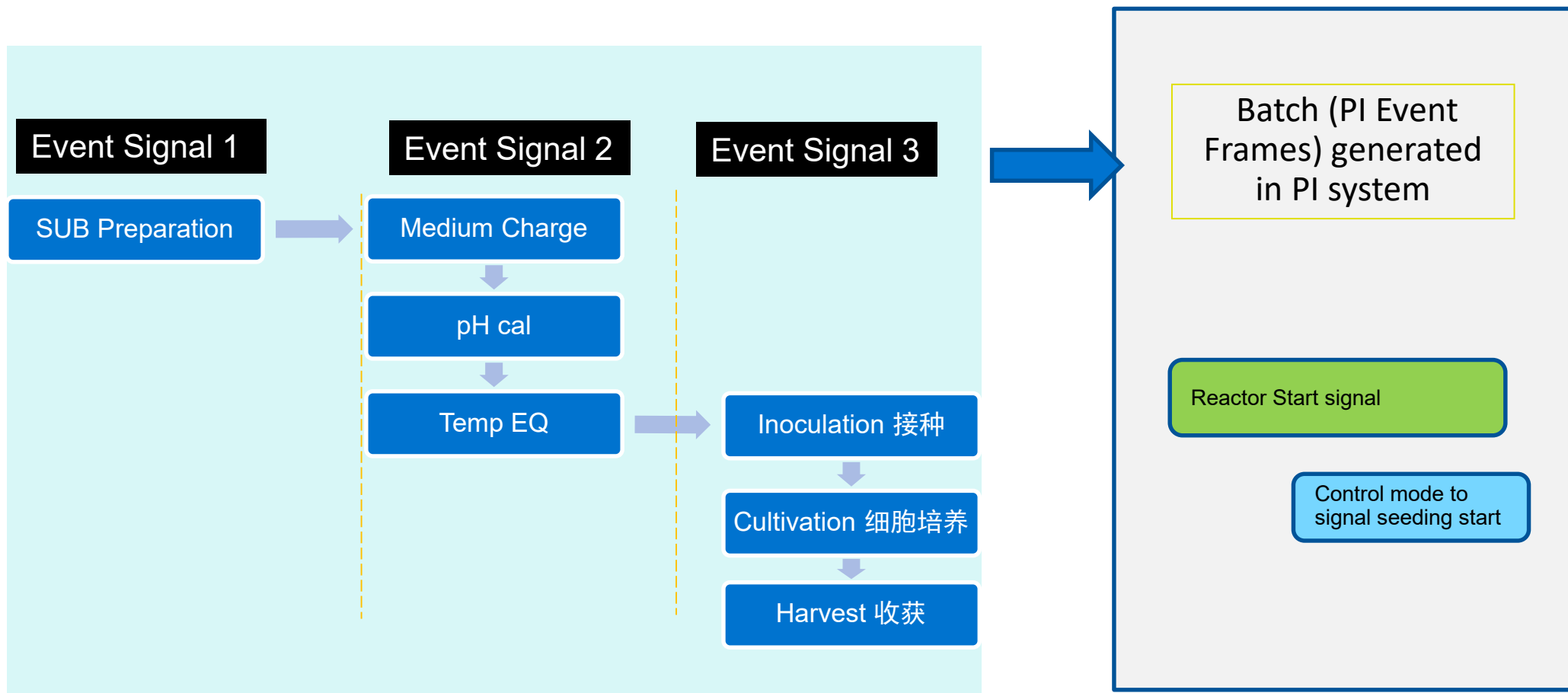


# 生物制药工艺: 本质上基于批次过程

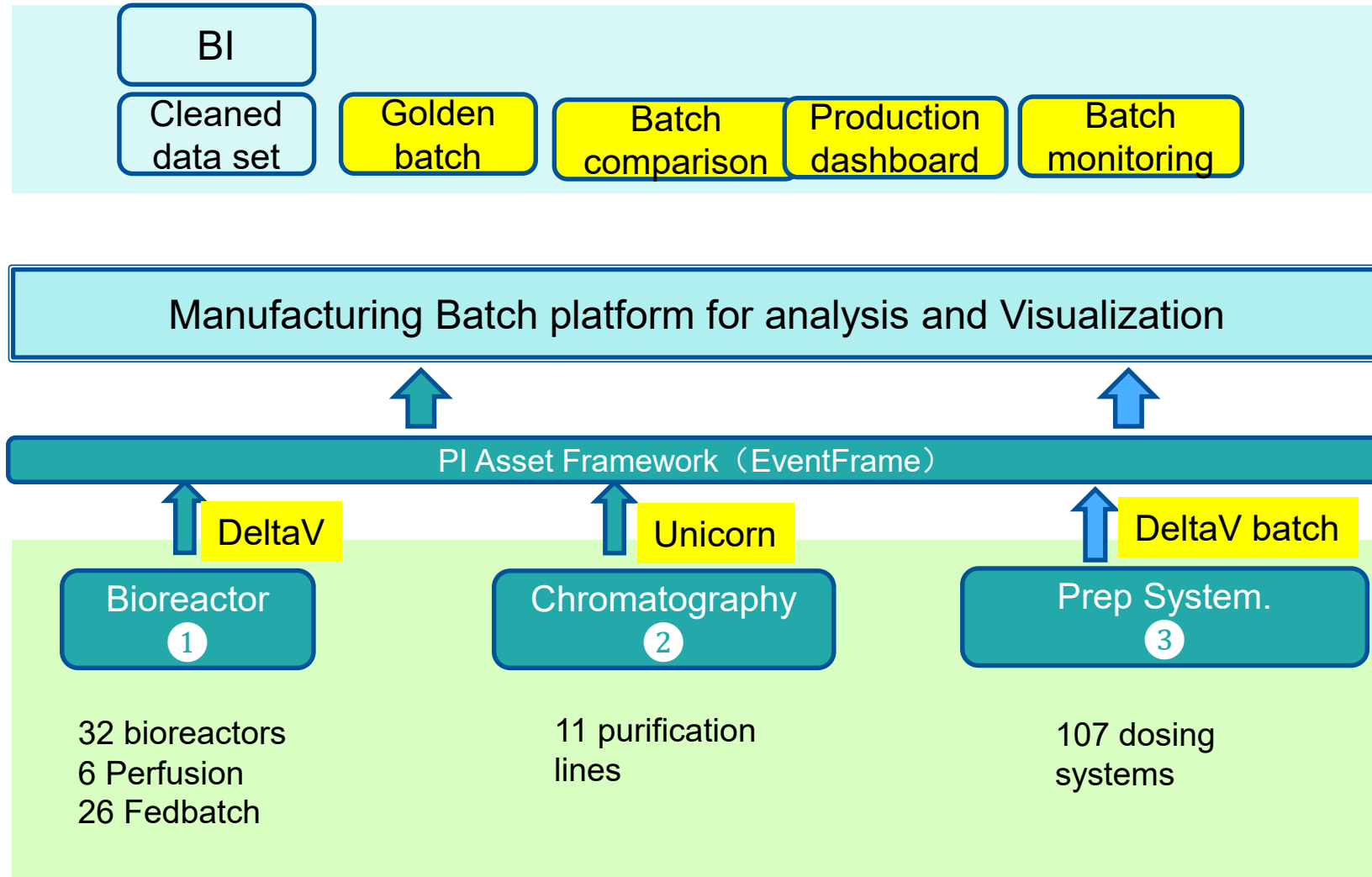
## Biopharmaceutical process: essentially based on a batch process



# 细胞培养流程的批次化 - Batch cell culture process

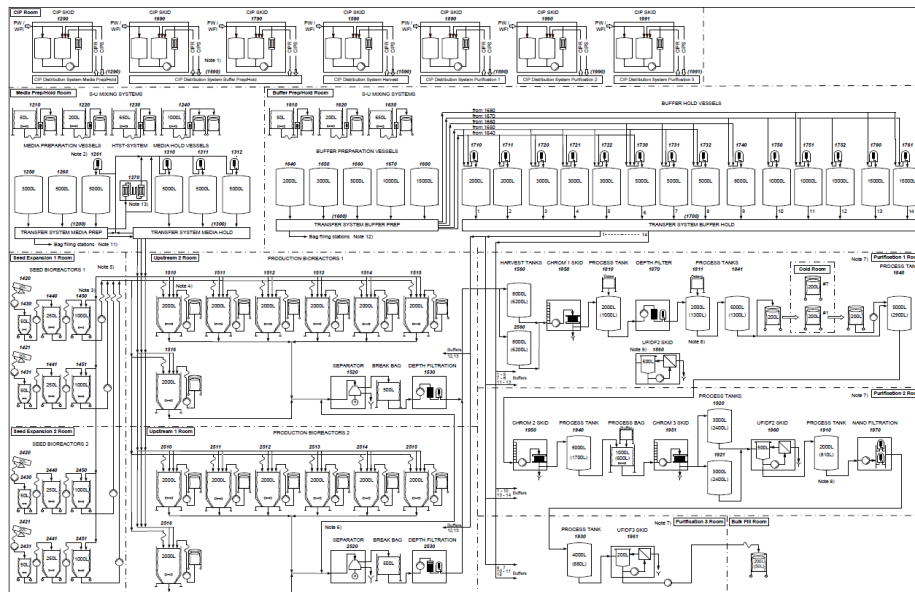


# 生物制药批次信息数据平台 - Biopharmaceutical batch information data platform



# SMART MAINTENANCE

- **Traditional: Preventive Maintenance (PM) & Repair**
  - PM as defined by GMP SOP, impacting production slot & cost
  - Repair when there is malfunctioning, leading to product loss
- **Smart Maintenance: Predictive Maintenance based on utilization and status**
  - Early prediction/detection of component failure
  - Reduce deviations caused by malfunctioning between PMs
  - Reduce loss of production slots and cost for maintenance



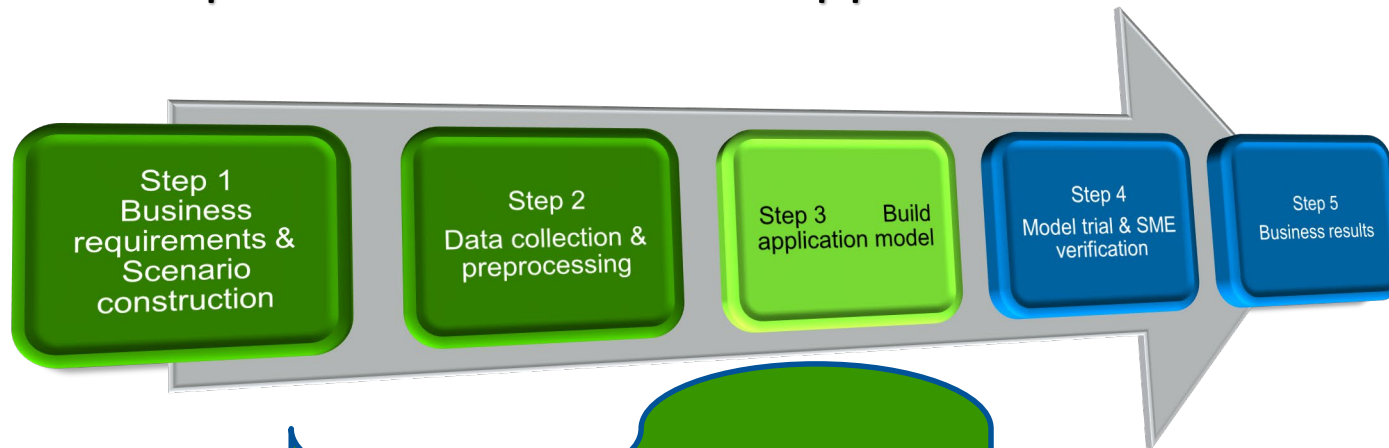
## Super Skids for Biologics Manufacturing

- **Vessel Systems**
- **CIP/ SIP Operations**
  - >200 CIP/SIP operations
  - >6000 automatic valves
  - >4000 gasket connections
- **Auto valves**
- **Gasket connections**
- **PM at fixed time interval**



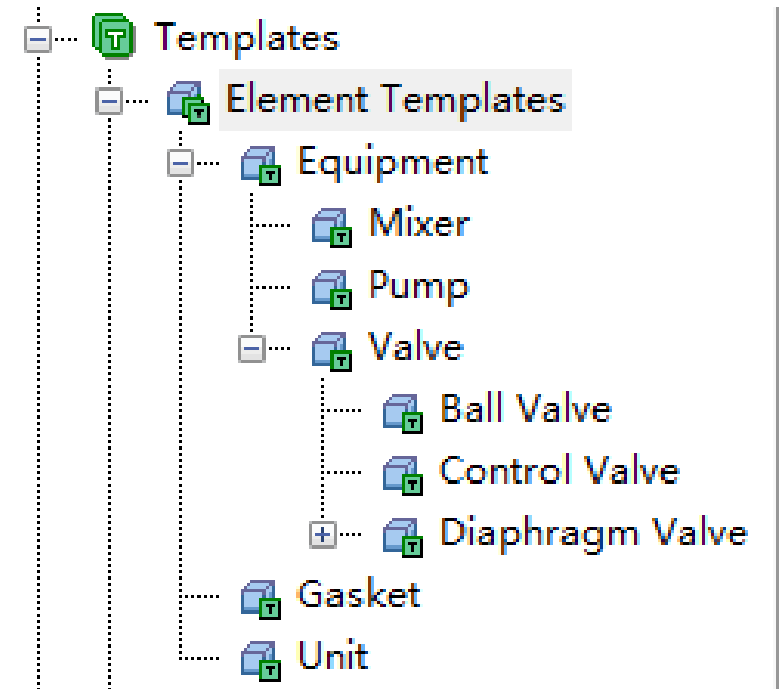
# Smart Maintenance System

## Five steps to build smart data application



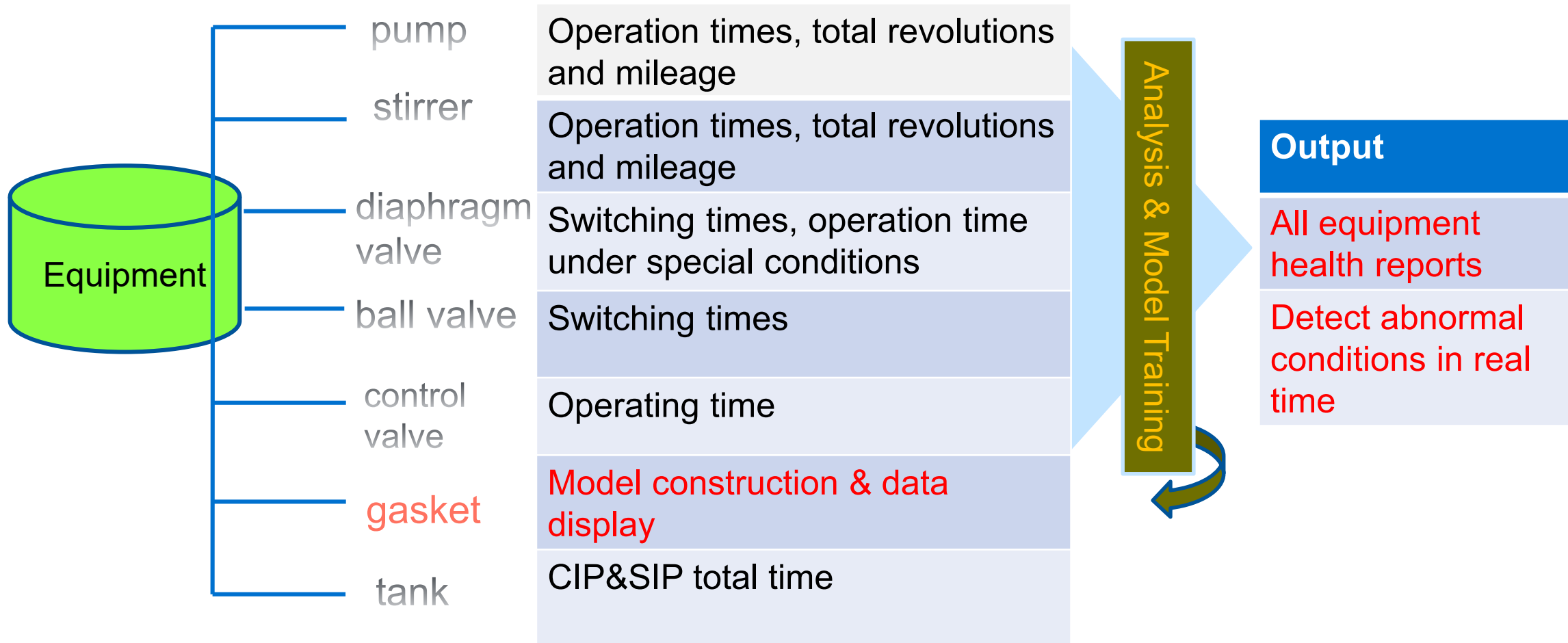
- PI System as the infrastructure
- PI AF is used to structure the plant's equipment assets
- The usage of each equipment component can be tracked according to different operating conditions

## Equipment model built by PI AF



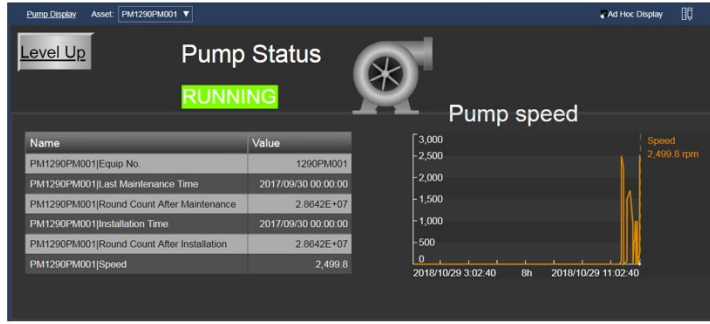
# Smart Maintenance System

Based on the operating status of the equipment in the operating cycle



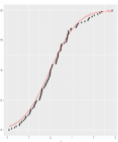
# Smart Maintenance Dashboard

- When was it installed?
- When was its last maintenance?
- Round Count – “the mileage”
- Operation counts – “Stop/go”
- Speed profile



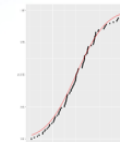
- ❖ Alert is generated when CHI exceeds set limit, calling for maintenance
- ❖ Alert is listed up based on its criticality
- ❖ Alert disappears when maintenance is performed
- ❖ The system now is in test bed for real world data

- Operation Count
- Round Count – “mileage”

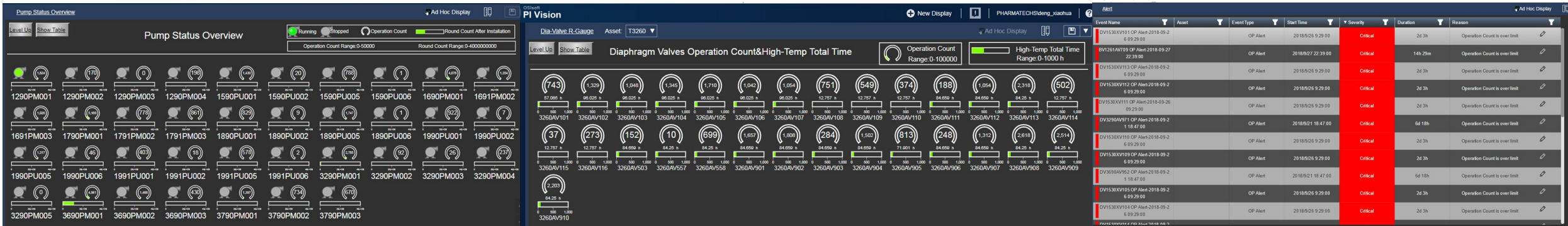


Aggregated Component Health Index (CHI)

- Operation Count
- High Temp Exposure Time



Aggregated Component Health Index (CHI)





# Achieved

- Batch process data platform to realize real-time data analysis and deviation warning
- Smart maintenance system, monitor the health coefficient of key components of liquid dispensing system in real time and guide equipment maintenance based on early warning
- Asset structures and data models can be seamlessly applied to global factories

## Next step...

- Aiming at the current smart maintenance system, to introduce AI tools (machine learning, model training) such as SIMCA, R and MATLAB
- Model trial and data verification to optimize the maintenance process
- Promote Wuxi's PI System to global production bases

# The largest user community for operations data in Life Sciences?

850+ members, 250+ organizations

[www.pisquare.osisoft.com/groups/life-sciences](http://www.pisquare.osisoft.com/groups/life-sciences)



## Steering Committee

- Biogen
- Lilly
- Genentech
- Roche
- Regeneron
- BioMarin
- DRL
- Novo Nordisk



# Faster and More Reliable Operations: A FUJIFILM Diosynth Biotechnologies' Case Study

## Challenge

Being able to review and verify process performance is a key step in biopharmaceutical manufacturing. However, traditional chromatography review, like visually reviewing elution peaks, was a **time-consuming, error prone, and paper-intensive** process.

## Solution

Paring Sartorius Data Analytics tools with OSIsoft PI technology FUJIFILM Diosynth Biotechnologies developed a new digital chromatogram dashboard for **enhanced chromatography review.**

## Benefits

**Shorter** review times, resource expenditure **optimized** tenfold, paper footprint **reduced** by ~10,000 sheets/year, **on demand data** accessibility, and **increased partner trust** and collaboration opportunities.

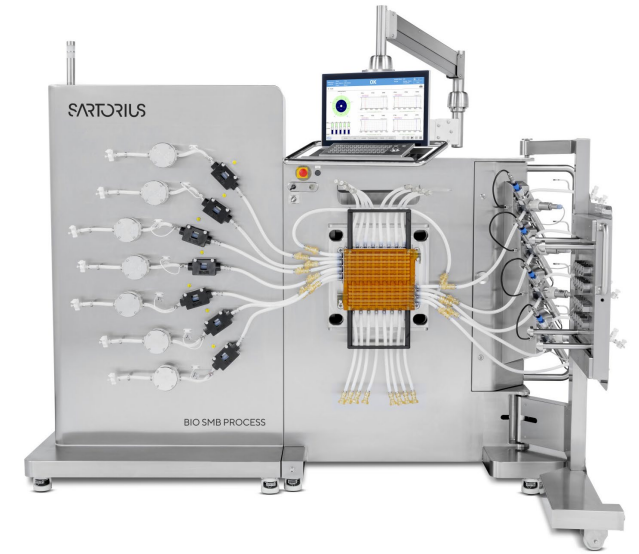
Do not duplicate or distribute without written permission from FUJIFILM Diosynth Biotechnologies.

# Using Data Analytics to Overcome Chromatography Review Challenges

Traditional chromatography process monitoring, like visually reviewing elution peaks, is time consuming, error-prone, and not fit for continuous bioprocessing.

Multivariate data analytics tools like SIMCA® and SIMCA®-online enhance chromatography monitoring and review by detecting small deviations in peak shapes before traditional methods can.

- Monitor process consistency
- Identify process failures and other trends before they become problematic
- Easily detect and correct column, cycle, or batch deviations
- Expedite chromatography release process - bringing you one step closer to RTRT

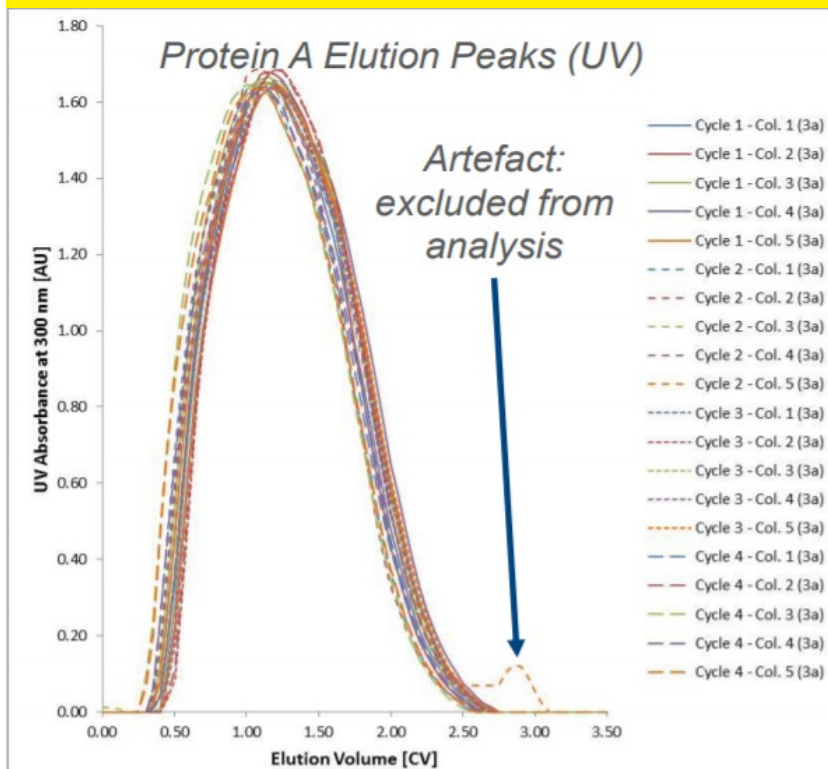


BIOSMB Process 80/350

Single-Use Continuous Chromatography Systems For Perfusion and Batch Bioreactor-based Processes

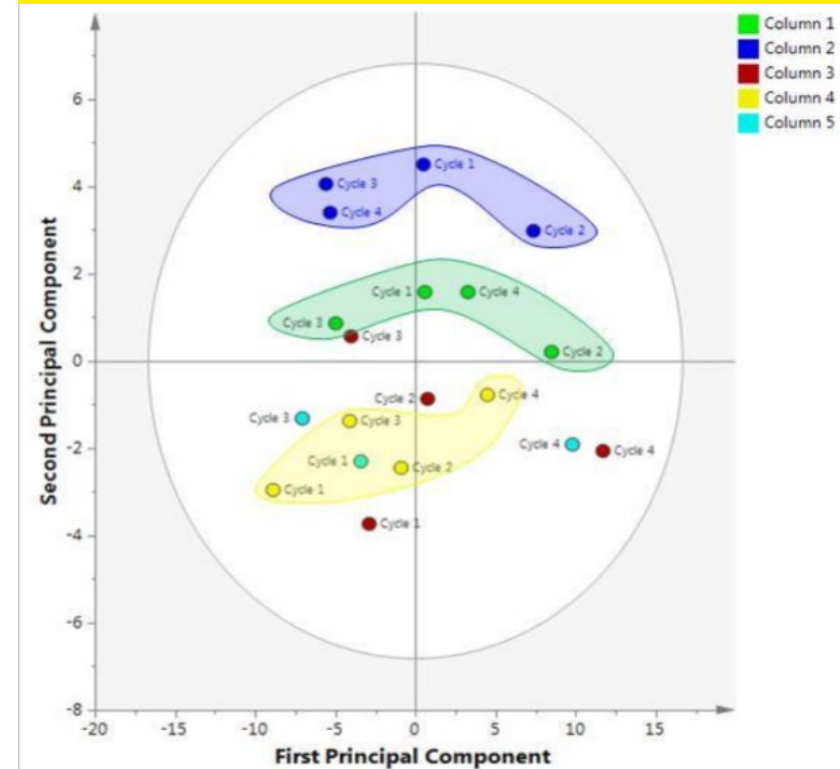
# Identify Column-to-Column Variations

## Traditional Univariate Analysis



Significant Variations Difficult to Detect

## Multivariate Analysis

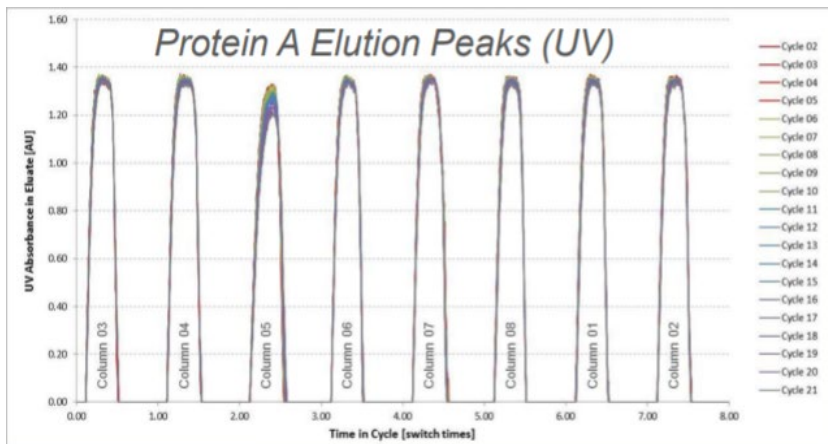
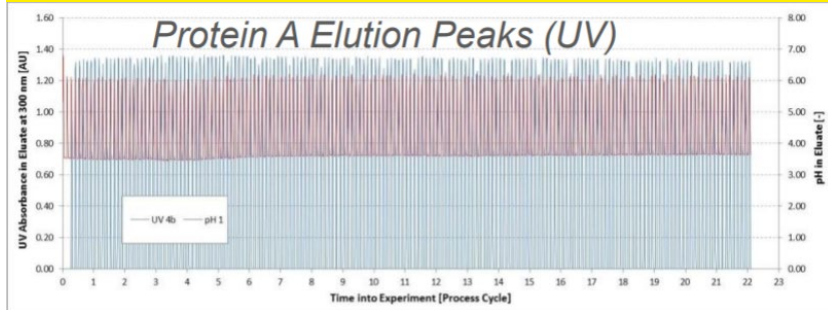


Column-to-column variations Easy to Detect



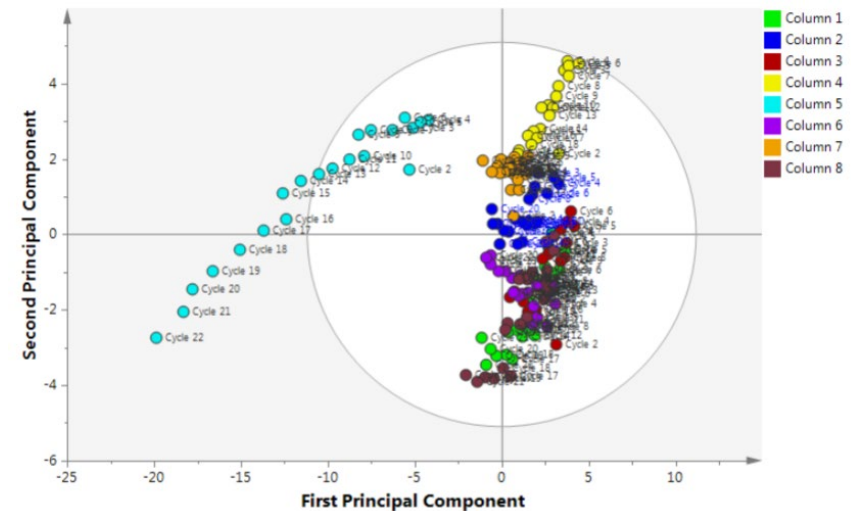
# Identify Column Malfunctioning

## Traditional Univariate Analysis



Cycle-to-cycle overlay shows some effect in column 5

## Multivariate Analysis

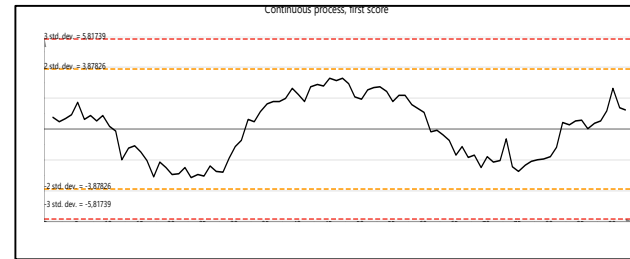


Performance decay in column 5 clearly indicate. Investigation revealed inadequate cleaning conditions.

# Implementing Model Based Batch Review



**1. Model Building:** Trained Model Builder Collects Historical Data (IPC, Analytical, Raw Material) and Builds Model with SIMCA®

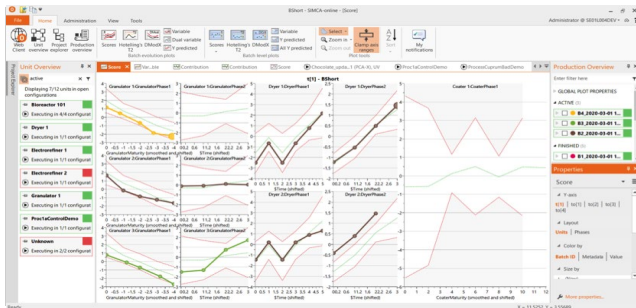


$$t[1] = x1*Temperature + x2*Pressure + x3*Speed + x4*pH + \dots$$

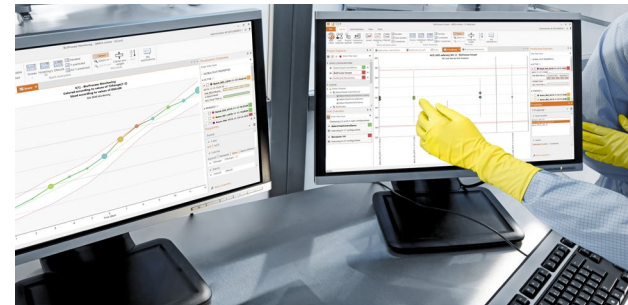
**2. MVDA Compression:** Large datasets compressed into smaller latent variables allows for a simpler representation



**3. Connecting Data Sources:** In order to enable real-time monitoring the model is connected to the original data sources



**4. Monitor Incoming Batches:** SIMCA®-online applies MVDA in real-time for monitoring new batches

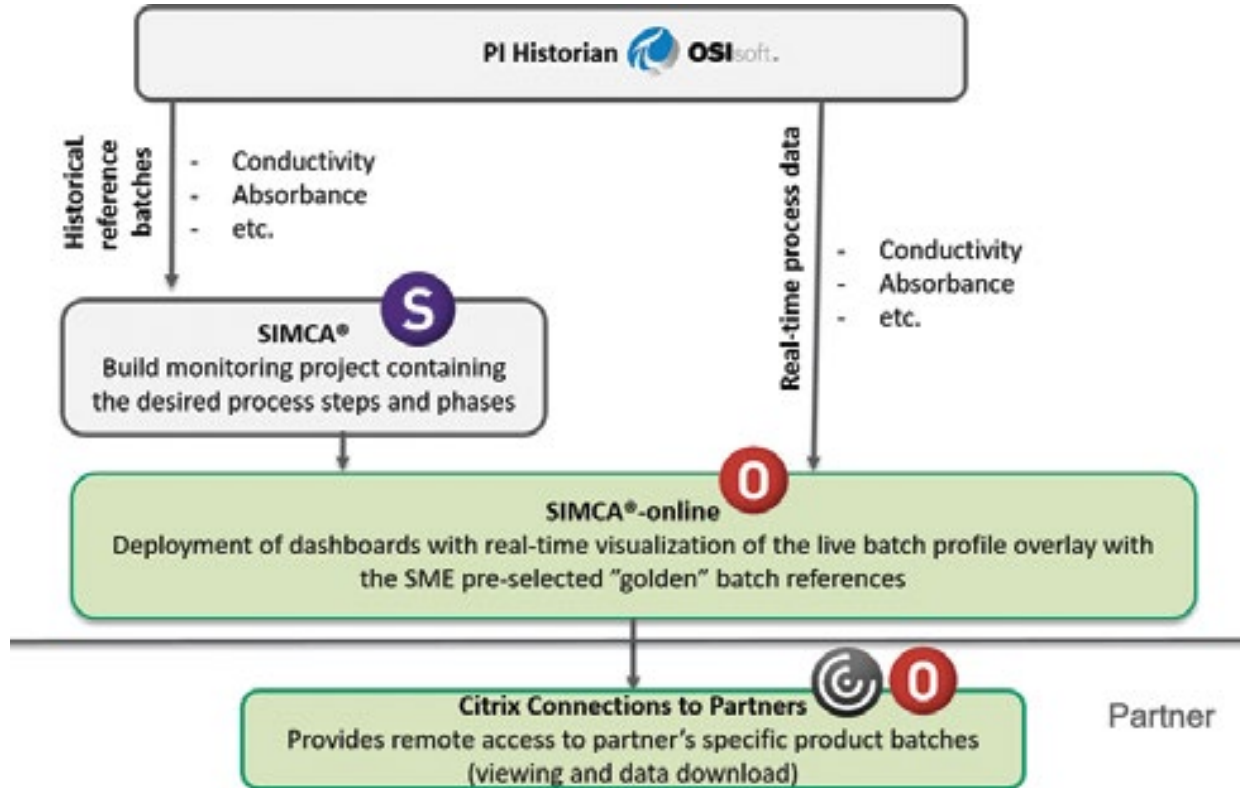


**5. Real-time Quality Assurance:** Trends are viewed live on plant floor or from control room



**6. Immediate Risk Mitigation:** alarms highlight process deviations and allowing for fault faster reaction

# System Set-up: A FUJIFILM Diosynth Biotechnologies' Case Study

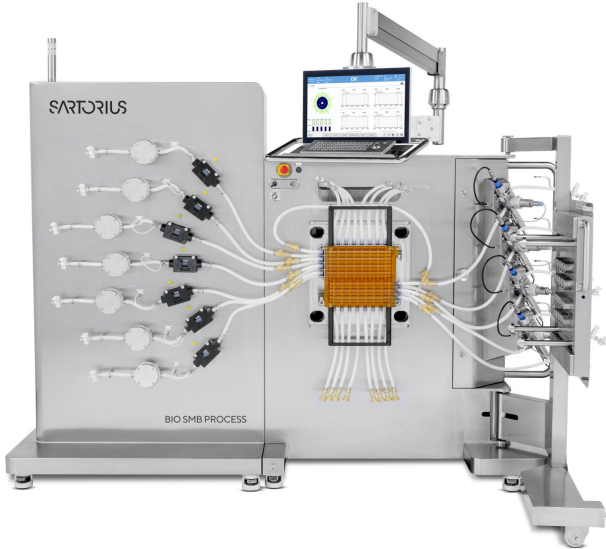


- Historical data stored in the OSI PI system were used as reference batches and to build the digital chromatogram project
- During production, data ingested from the manufacturing SCADA system to OSI PI system were transmitted to SIMCA®-online through a SimAPI at real-time frequency
- The visualization dashboard were then made available to operators, quality assurance personnel reviews, and the partner

Do not duplicate or distribute without written permission from FUJIFILM Diosynth Biotechnologies.

# Improving Partner Data Visibility

CDMO Process



Real-time Process Window and Data Download



Partner



# Benefits Realized: A FUJIFILM Diosynth Biotechnologies' Case Study

- **Shorter** review times

Accelerated Delivery

- Resource expenditure **optimized** tenfold

COGs Optimization

- Paper footprint **reduced** by ~10,000 sheets/year

Positive Environmental Impact

- **On demand** data accessibility

Improved Process Traceability

- **Increased** partner trust and collaboration opportunities

Increased Reliability

Do not duplicate or distribute without written permission from FUJIFILM Diosynth Biotechnologies.

# CDMOs are Putting a Greater Importance on Data and Data Analytics

“Data integrity deficiencies are cited in **65% of all FDA warning letters**”<sup>1</sup>

Consequently this has put pressure on CDMOs to implement data management systems that provide **transparency, preserve customer confidentiality**, and meet **regulatory guidelines**

CDMOs have realized the **additional value opportunity** that comes with adopting data management systems along with data analytics tools and have transformed these new technologies into **service offerings**



SARTORIUS

# Acknowledgements and References

Thanks to **Martin D. Jensen** and **Ricardo F. Caroço** from FUJIFILM Diosynth Biotechnologies in Hillerød, Denmark

Enabling Digital Chromatogram Review for a Faster and More Reliable Operation, M. Jensen, R. Caroço, *Bioprocess International*, Industry Innovators, pg. 30, 2020-2021



FUJIFILM Diosynth Biotechnologies

**DOWNSTREAM PROCESSING Chromatography**

## Enabling Digital Chromatogram Review for a Faster and More Reliable Operation

**FUJIFILM**  
Diosynth biotechnologies

Chromatogram review is a monitoring method used to verify process performance in packed-bed chromatography processes. By observing key process parameters such as chromatography column outlet conductivity or UV absorbance, it is possible to identify the signs of a poorly packed column, resin degradation, or equipment malfunction. Therefore, chromatogram review is implemented as an in-process control (IPC) to decrease variability and identify suboptimal performance, thereby enhancing yield and ensuring high product quality (1).

The industry standard practice relies on trending univariate parameters (e.g., chromatographic peak asymmetry and product yields) and on performing a qualitative visual comparison of chromatography profiles against a reference batch. The latter is set as a control check to be performed by operators before proceeding with the next stage of a process.

That assessment used to be executed by printing a physical copy of the chromatogram of a batch and then comparing that printout with a standard operating procedure (SOP) that contains the reference chromatogram. The same process would be reviewed later by quality assurance personnel. That approach led to a time-consuming, paper-intensive process that was prone to mistakes caused by the variability inherent to visual comparison of chromatogram profiles not only in different plots, but also on different sheets of paper.

Now that process has been reimagined as part of the drive toward Industry 4.0 and the desire to become a digital facility, enabled with systems that can respond to changes in real-time and act proactively with the necessary corrective behaviors.

With the new digital chromatogram dashboard, printing becomes unnecessary, and multiple phases and parameters are viewed simultaneously, leading to shorter review time. Reliability is increased because the dashboard enables monitoring of a live batch overlaid on reference batches preselected by a subject matter expert. That is facilitated by implementing the SIMCA® and SIMCA®-online software suites by Sartorius (2). As a leading contract development and manufacturing organization (CDMO), Fujifilm Diosynth Biotechnologies strives for increased partner trust and collaboration. Thus, these dashboards (and many others with multivariate models) can be accessed by a partner, providing it with a real-time window to the process (Figure 1).

By digitalizing the existent business process and

Figure 1: System and data communication diagram: historical data available in the data historian are used as reference batches and to build the digital chromatogram project. During production, data ingested from the manufacturing SCADA system to the data historian (DS) are transmitted to SIMCA-Online through SimApi at real-time frequency (2). The visualization dashboards are made available to operators, quality assurance personnel reviewers, and the partner.

expenditure is optimized tenfold (shorter review time, fewer investigations expected), the paper footprint is reduced (~10,000 sheets/year), and data accessibility is on demand. The review process is not hidden and is readily available to different departments.

**REFERENCES**

- 1 ICH Q7: Good Manufacturing Practice Guidance for Active Pharmaceutical Ingredients. The International Council for Harmonisation of Technical Requirements for Pharmaceuticals for Human Use (ICH), 2016.
- 2 Data Analytics AB. SIMCA-Online Technical Guide. Sartorius Stedim, 2020; <https://umetrics.com/kb/simca-online-technical-guide>.

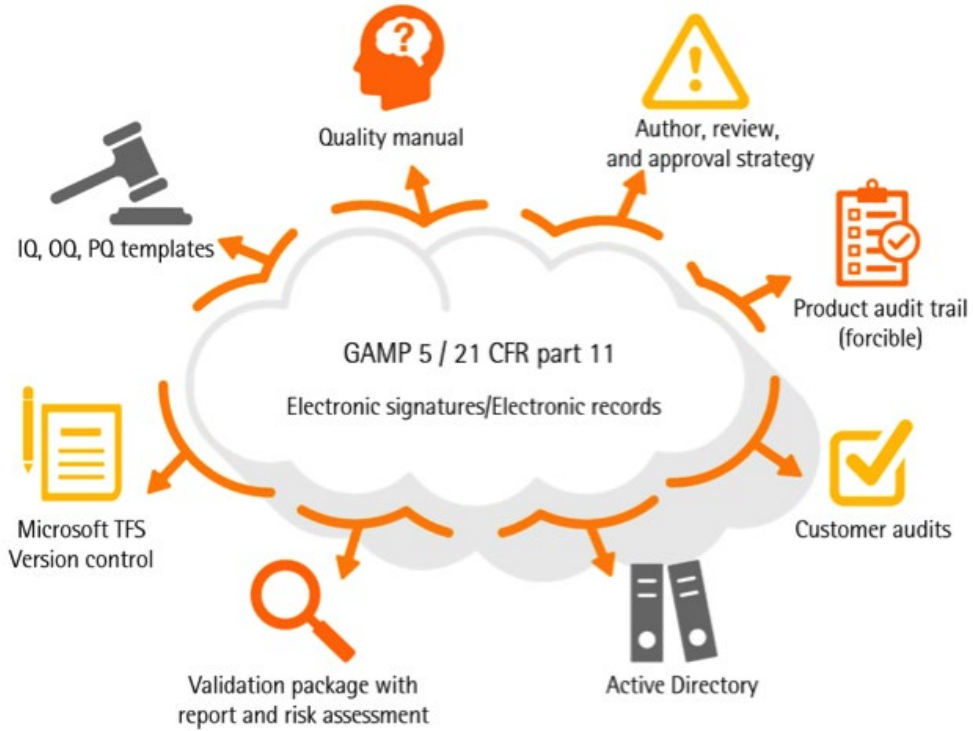
**Martin D. Jensen** is engineer III, manufacturing services, and **Ricardo F. Caroço** is a process analytics engineer at Fujifilm Diosynth Biotechnologies.

Enabling Digital Chromatogram Review and More Reliable Operation

# Interested in Reading More about Data Integrity?

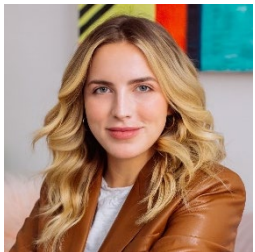


[How to ensure data integrity and compliance of your data analytics systems](#), Erik Renberg (2018)





# Thank You!



Tiffany McLeod, Market  
Manager Pharma  
Biopharma,  
Sartorius Data Analytics



Petter Mörée,  
Industry Principal,  
Global Lifesciences,  
OSIsoft

**SARTORIUS**