



Capturing the value of Continuous Bioprocessing through MVDA

*Marc Bisschops
Principal Scientist
Pall Life Sciences*

This presentation is the confidential work product of Pall Corporation and no portion of this presentation may be copied, published, performed, or redistributed without the express written authority of a Pall corporate officer

© 2015 Pall Corporation

Umetrics User Group Meeting
Princeton (NJ), September 9, 2015



Biopharmaceutical Manufacturing

- Biopharmaceutical industry is relatively young in terms of manufacturing operations
- Not all product attributes can be measured (at any given time)
 “The process is the product”
- Process validation is an inherent part of product quality control
- Dominated by batch manufacturing

... but changes are coming!

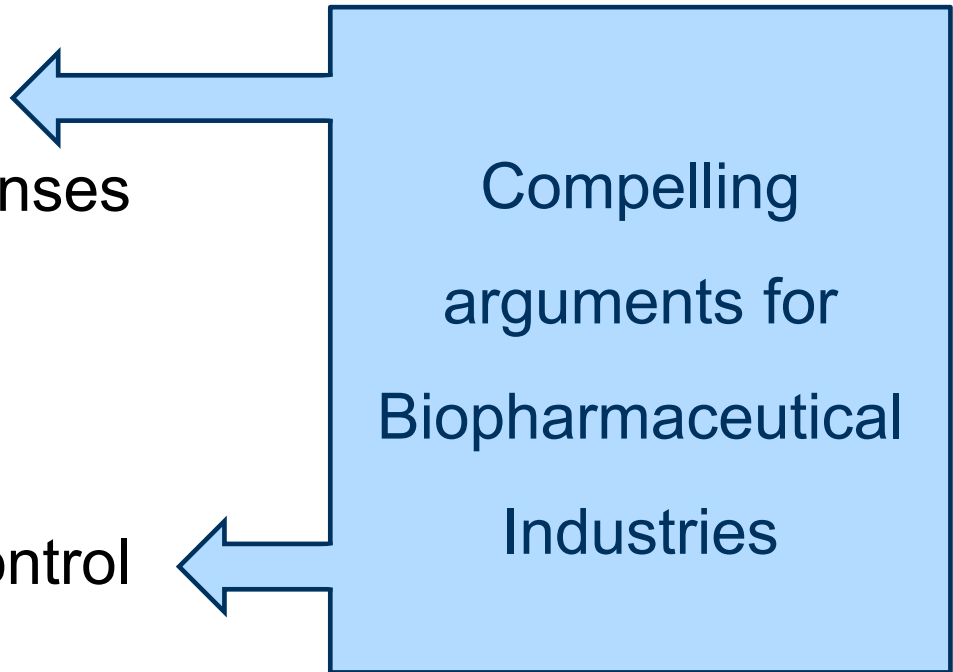


Continuous Manufacturing

Why are many mature industries moving towards continuous processing:

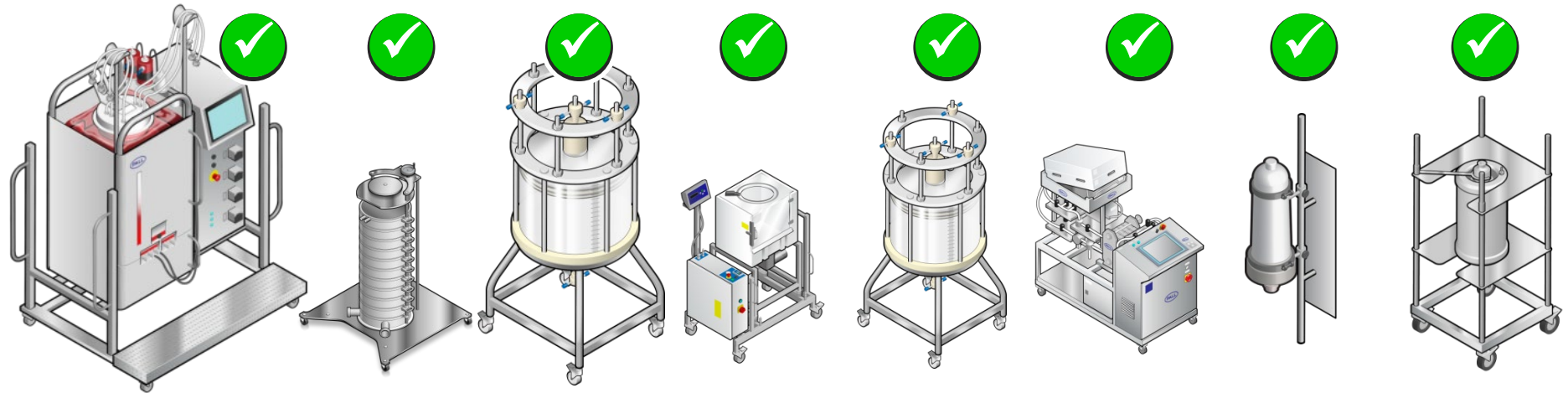
- Process economics:
 - improved capital utilization
 - reduction of operating expenses

- Process control:
 - Improved safety
 - Improved product quality control

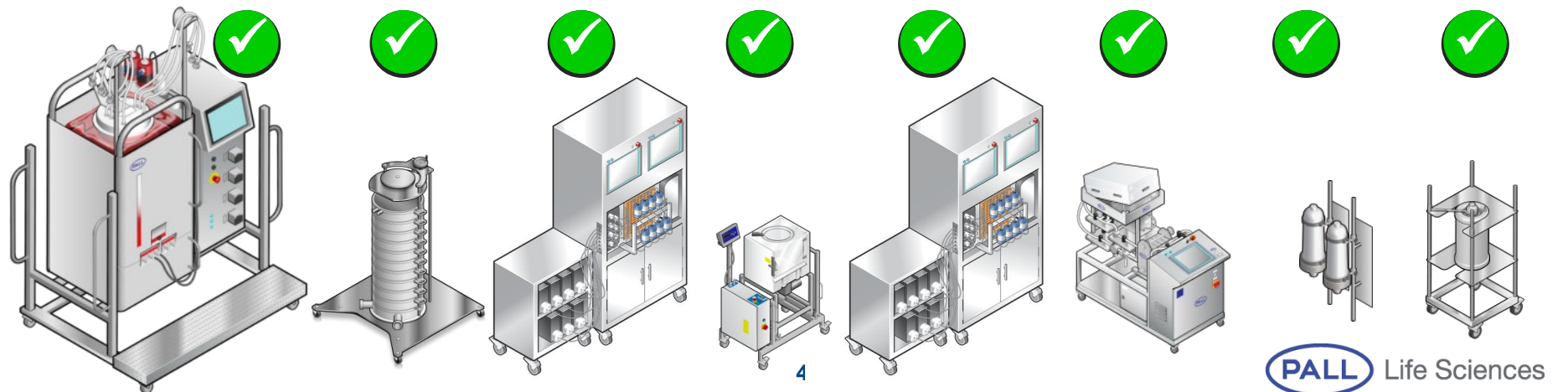


Integrated Continuous Bioprocessing

Fed Batch USP + Batch DSP

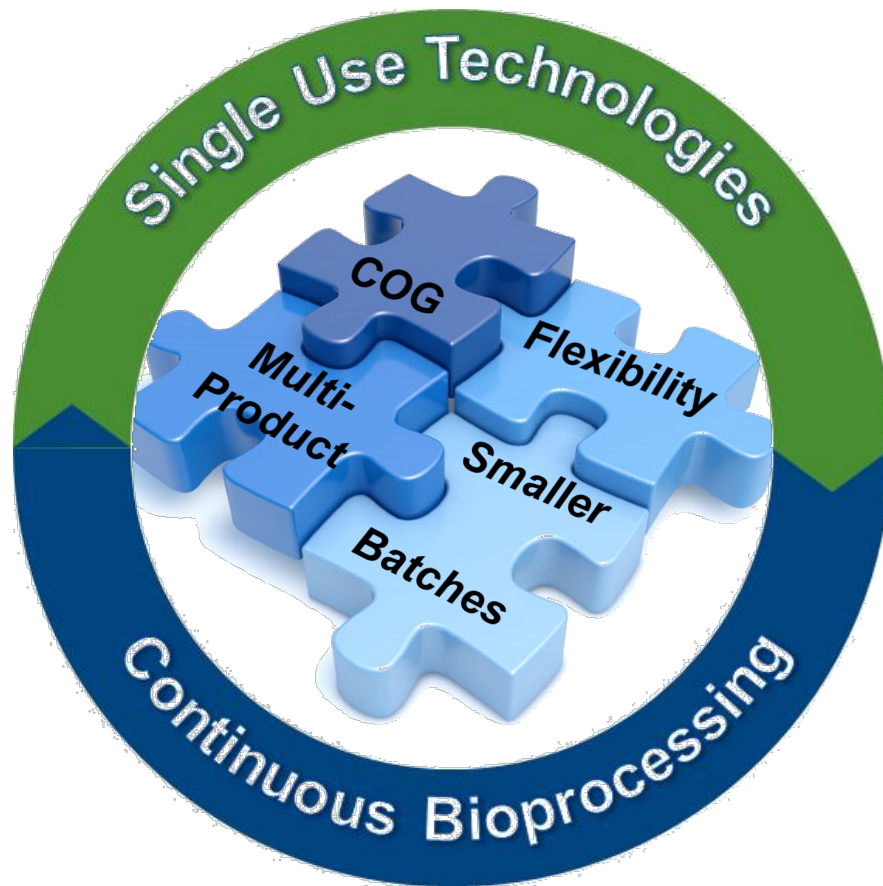


Fed Batch USP + Continuous DSP



Trends in Continuous Bioprocessing

Industry Trends



Considerations:

Most DSP steps are compatible with continuous bioprocessing:

- Filtration steps
- Flow-through chromatography steps

Key Enabling Technology:

- Multicolumn Chromatography (e.g. BioSMB[®]) for bind/elute chromatography

BioSMB® Technology

Key Benefits:

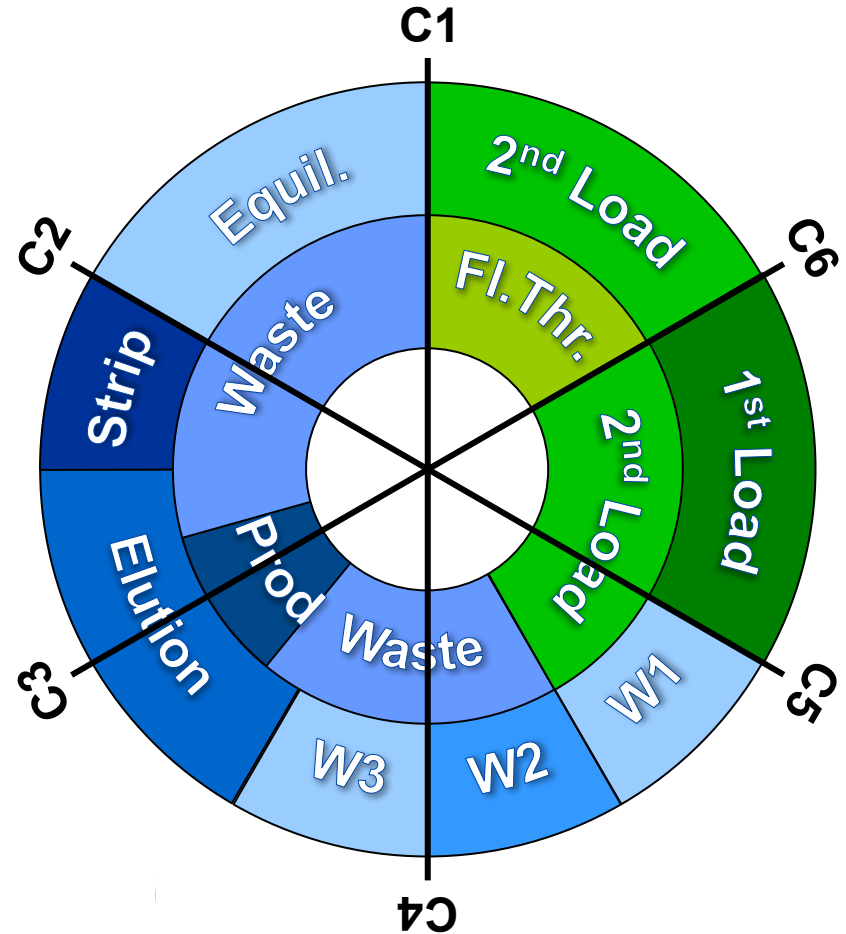
- Improved specific productivity
- Improved utilization of resin capacity
- Significant reduction in buffer consumption
- Enabler for integrated continuous bioprocessing
- Enabler for integrated single-use manufacturing



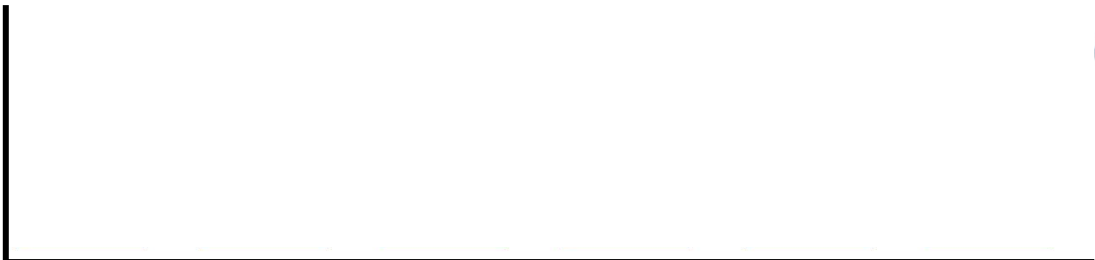
BioSMB[®] Technology

Highlights:

- Multiple columns work together to allow continuous feed
- Columns travel through the process (or actually vice versa)
- Each column results in one elution peak every cycle



UV Absorbance in Product Outlet





Evaluation of Chromatographic Performance

Traditional chromatography process monitoring:

Process performance monitoring	Comment
Column characterization (HETP and asymmetry)	Prior to process start
Critical parameters (pool volumes, yields, etc.)	Off-line analyses
Review of chromatographic peaks: <ul style="list-style-type: none">• Visual review• Moment analysis	Based on on-line data

**This Strategy may need to be reconsidered
for Continuous Bioprocessing**



Using old Tools for a new Approach

Multivariate Data Analysis (MVDA):

- A mathematical tool for data reduction
- Very strong for recognizing patterns in large and complex datasets

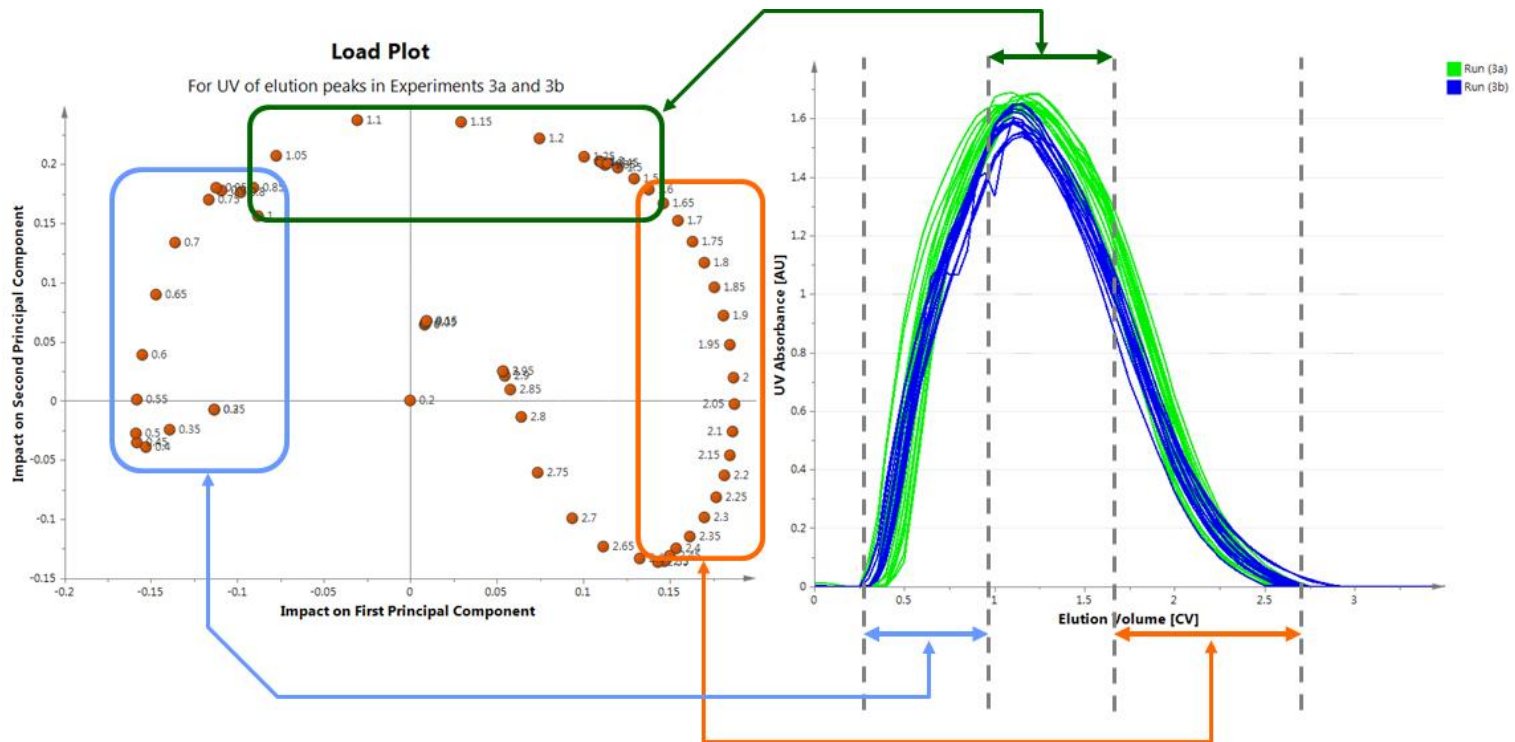
In this study, we will limit the MVDA to:

Principal Components Analysis (PCA):

- Not new: PCA was first described (invented) in 1901
- Mathematical transformation of data into orthogonal (principal) components

PCA for Chromatography

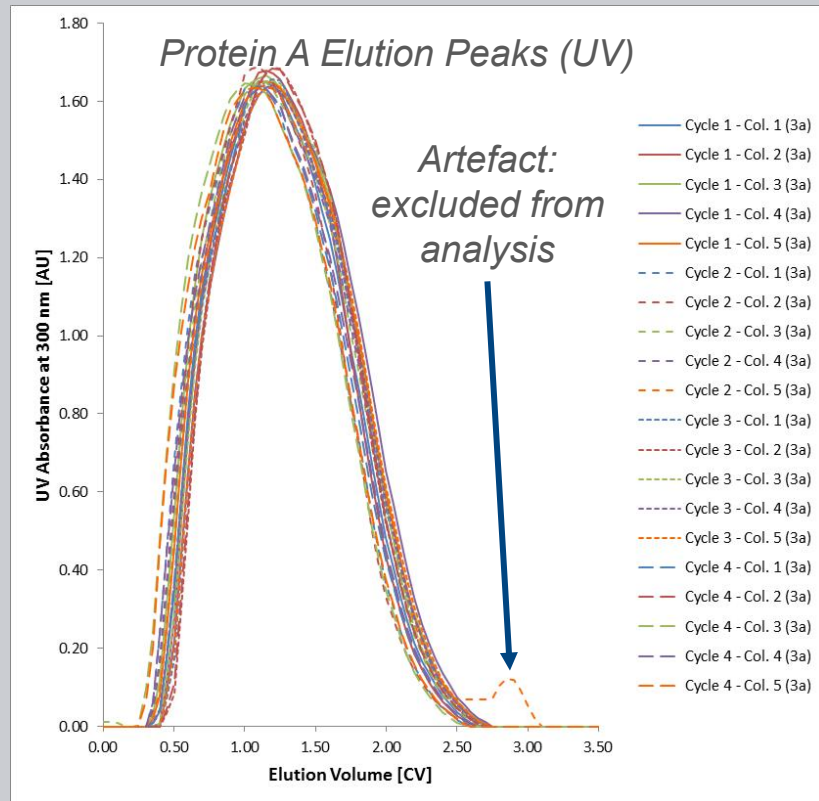
The Principal Components do not (necessarily) represent a physical characteristic of the chromatography peak, but...



The correlation between load plot and chromatography peak changes with the data set.

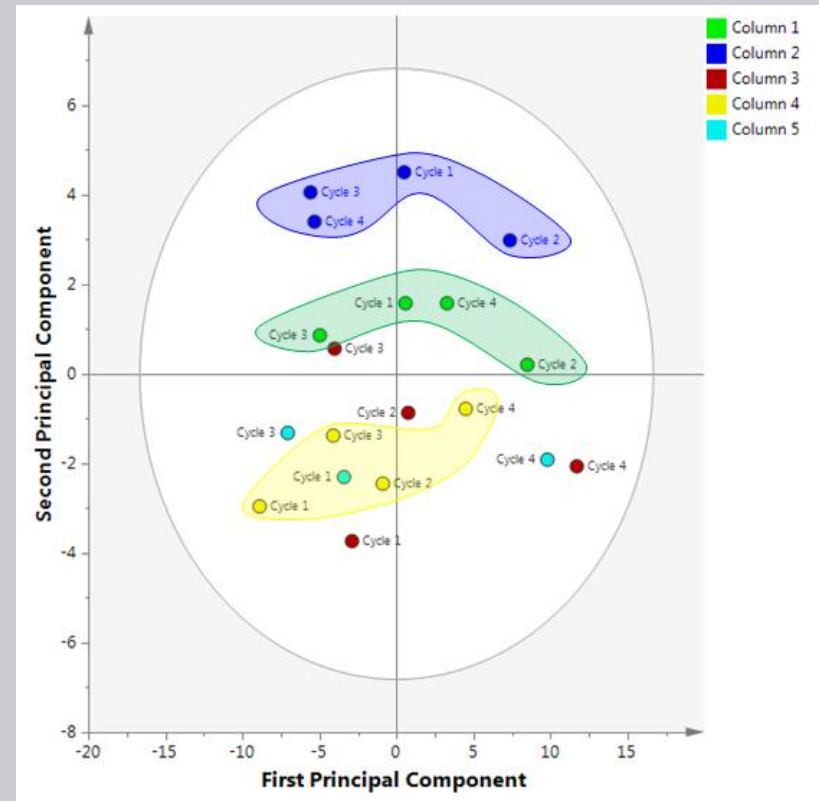
Case Study 1: Column to Column Variations

Monovariate Analysis



No significant variations detectable

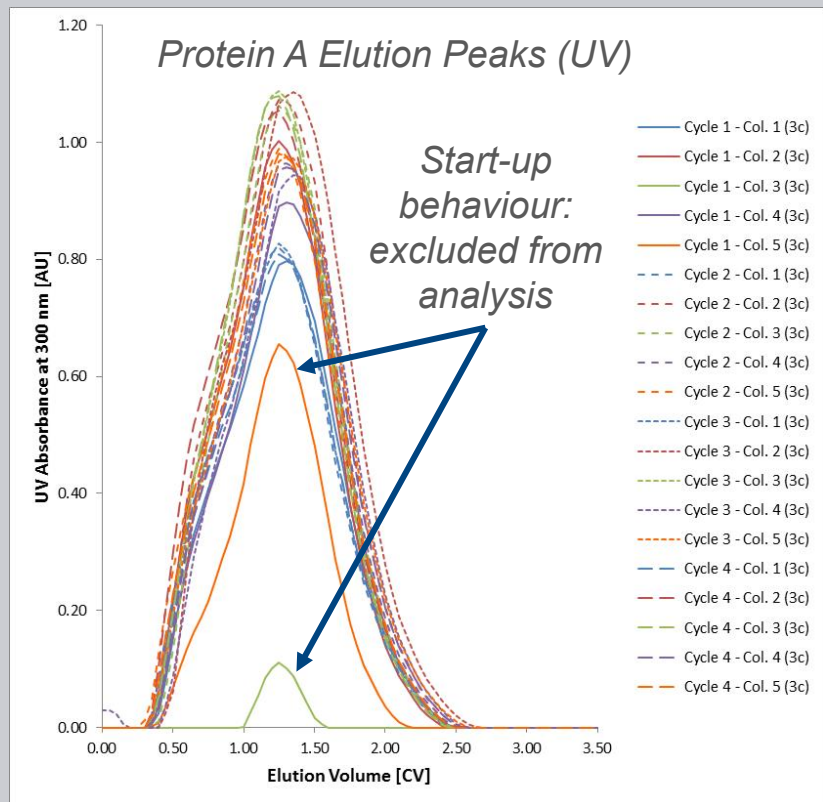
Multivariate Analysis



Main source for variation: Column-to-column variations

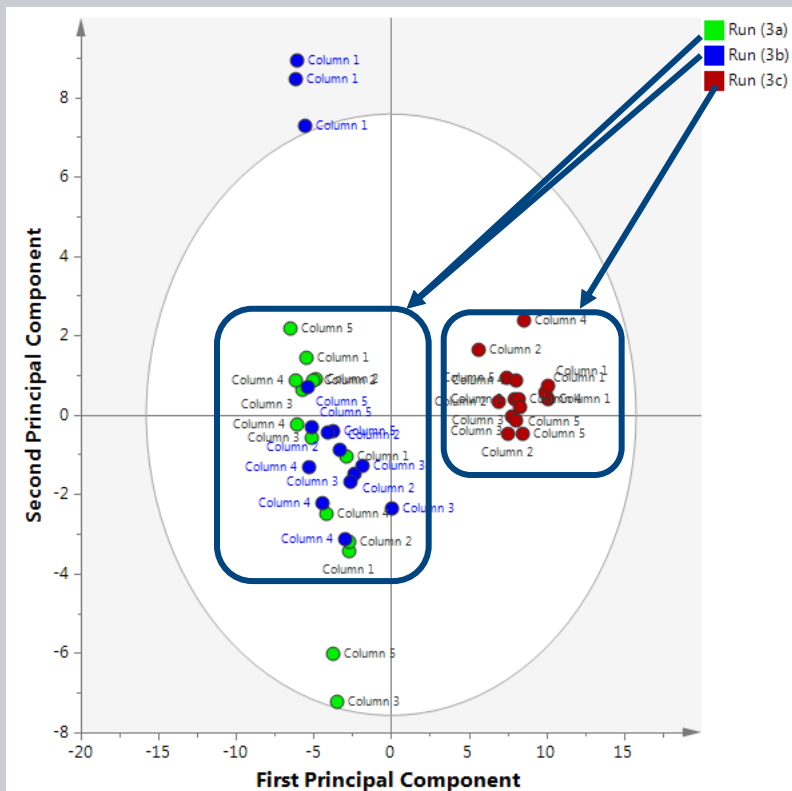
Case Study 1: Batch to Batch variations

Monovariate Analysis



No significant variations detectable

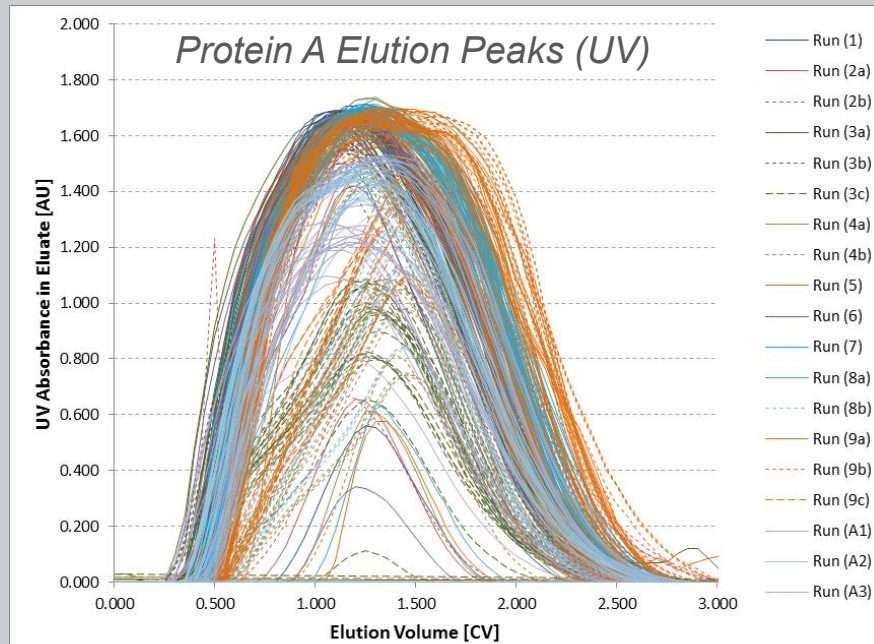
Multivariate Analysis



Reasonable consistency for two, but large variation for third batch

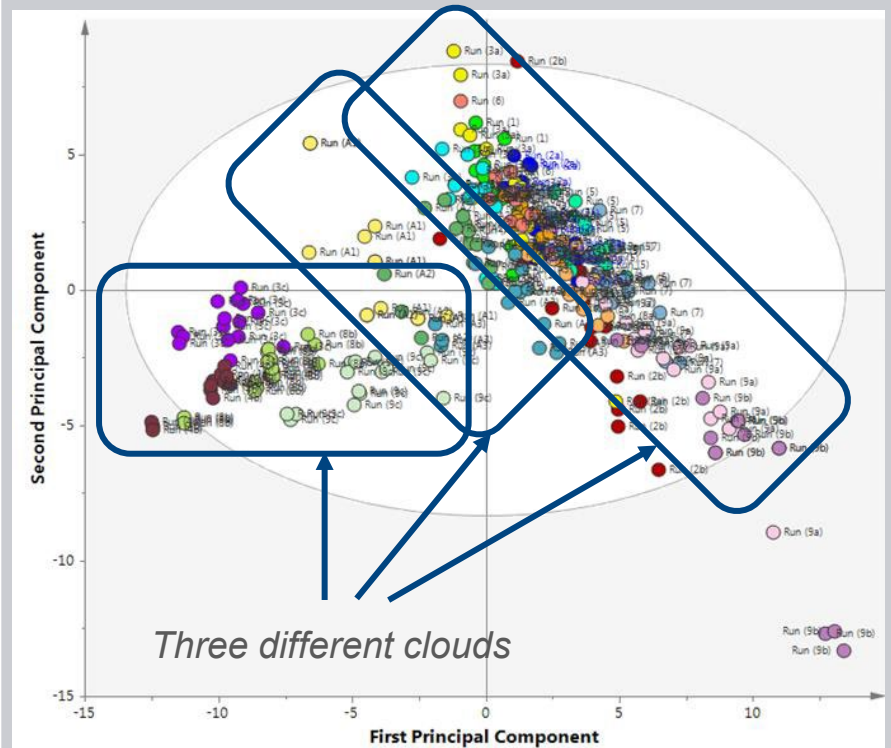
Case Study 1: Batch to Batch variations

Monovariate Analysis



Too much data to make any sense

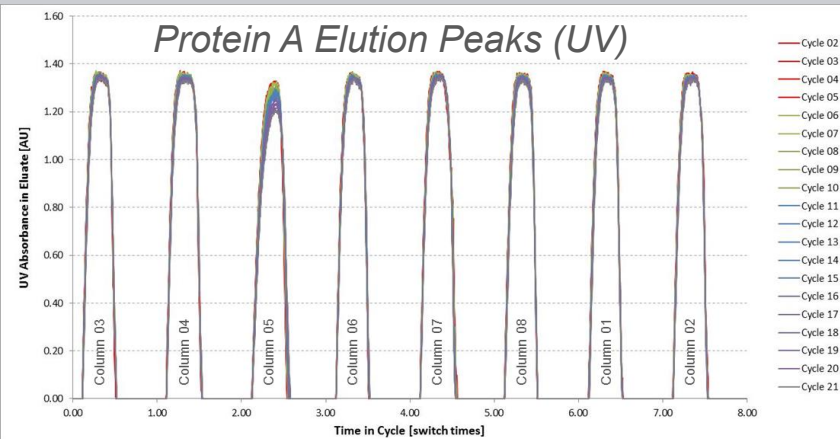
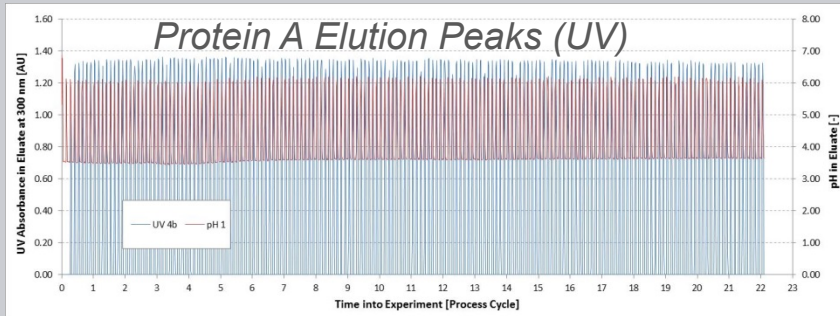
Multivariate Analysis



Harvest procedures impacted the PCA of PrA elution peaks

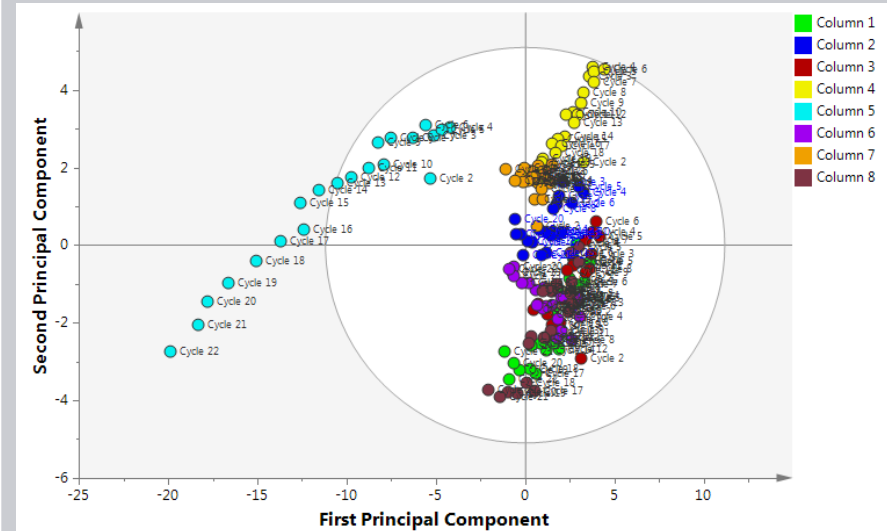
Case Study 2: Column Malfunctioning

Monivariate Analysis



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

Multivariate Analysis



Note: Performance decay in Column 5 was most likely related to inadequate cleaning conditions (not to the separation and/or technology itself)

Column 5 shows deviations from start of run



Opportunities

MVDA offers numerous opportunities for monitoring process consistency:

Opportunity (examples only)	Potential Approach
On-line column characterization	PCA on ΔP across columns (e.g. during equilibration or wash step)
Monitoring bed packing consistency	PCA on conductivity as column moves through different wash steps

Forward looking:

- Correlating response from MVDA to product attributes (QCA's) will bring us one step closer to parametric release
- Integrated process control (Process Analytical Technologies)



Conclusions

- Multivariate Data Analysis (MVDA) turns large datasets into (visual) information, thereby capturing the value of continuous bioprocessing
- Principal Components Analysis (PCA) can detect small deviations in peak shapes before traditional methods can:
 - Monitor process consistency (cycle-to-cycle reproducibility)
 - Detect column-to-column variations
 - Detect column failures and other trends before they becomes problematic

MVDA will help end-users providing evidence that they're in control of their process



Acknowledgements

Boehringer-Ingelheim (Fremont, CA):

- Jonathan Coffman
- Brian To
- Mark Strawn

Sanofi-Aventis (Frankfurt, GER):

- Annette Berg
- Hendrik Flato
- Stefan Schleper

Pall Life Sciences (Westborough, MA):

- Lynne Fick
- Steve Allen
- Ken Frank

