



System Performance Check (SPC): The Essential Partner for Multi-Attribute Method (MAM)

Jahziel Chase¹, Melissa Sato², Jared Auclair¹, Mike Knierman², Gordon Slysz²

¹Northeastern University, Boston, MA ²Agilent Technologies, Santa Clara, CA

Background and Hypothesis

Traditional quality control (QC) methods in biopharmaceutical analysis are time-consuming, resource-intensive, and often lack real-time insights. To address these limitations, a next-generation platform currently in beta testing is being developed to enhance Multi-Attribute Method (MAM) workflows. This platform integrates advanced LC-MS-based MAM with automated system performance checks (SPC), intelligent anomaly detection, and real-time monitoring of targets and critical quality attributes (CQAs). By improving specificity, compliance, and workflow efficiency, it aims to set a new standard for analytical rigor in biopharmaceutical QC.

We hypothesize that the integration of SPC and anomaly detection within MAM workflows will improve data quality, instrument reliability, and decision-making efficiency, enabling more robust and compliant biopharmaceutical characterization.

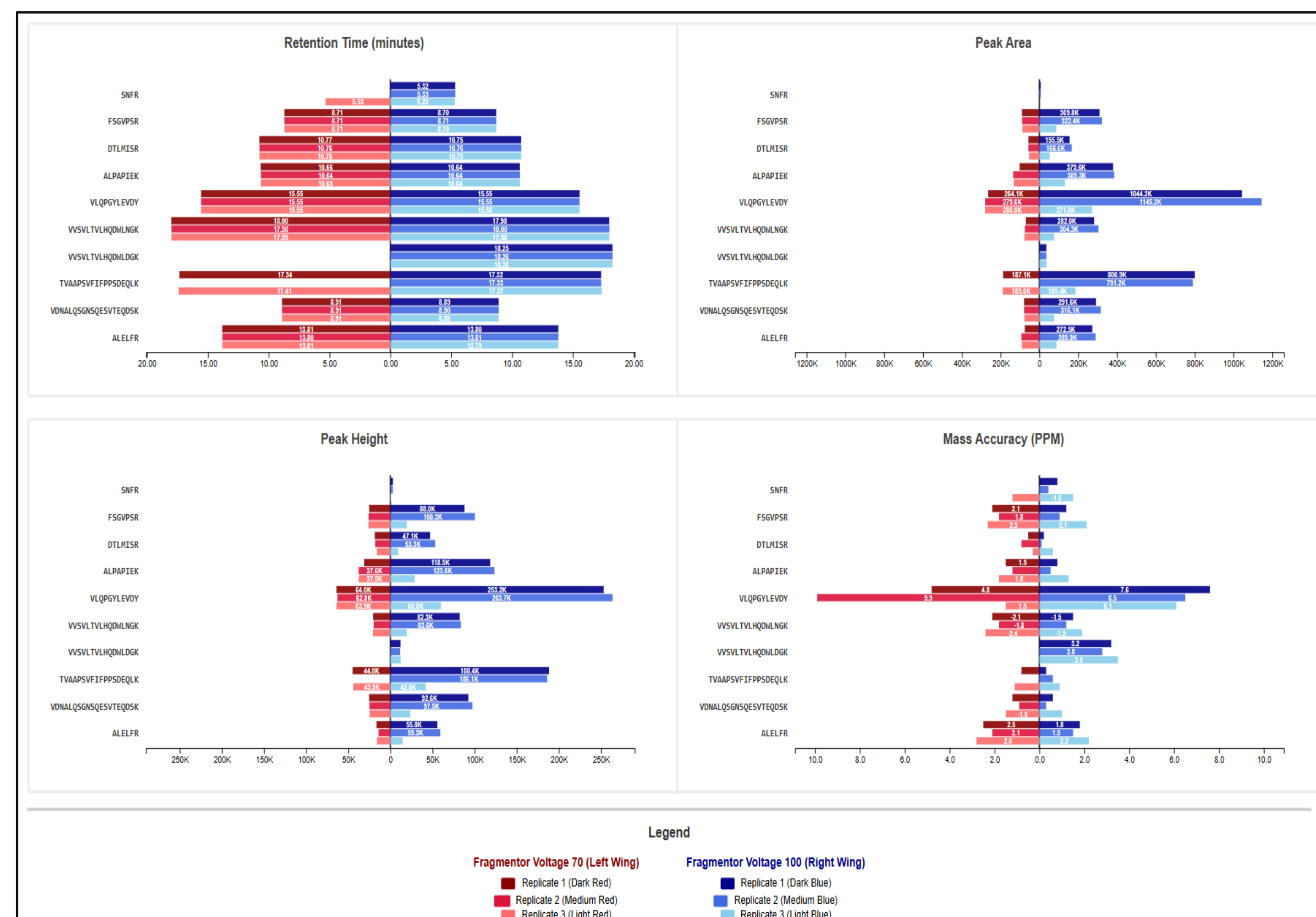
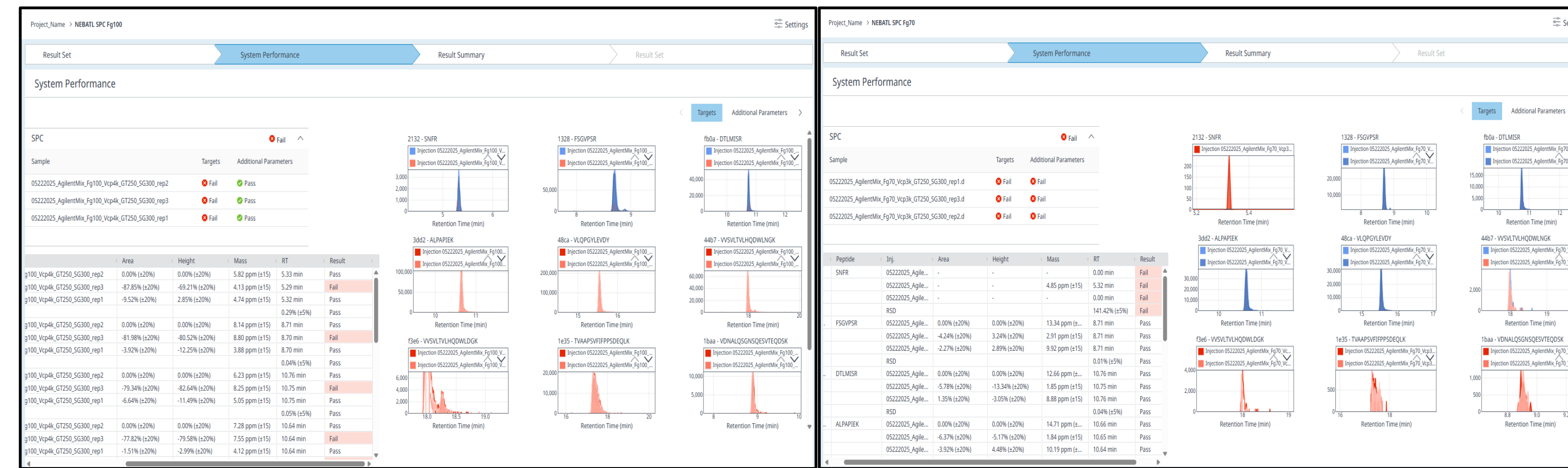
Objectives

- Identify limitations in conventional QC workflows.
- Demonstrate the role of SPC in ensuring reliable and reproducible MAM results.
- Evaluate the beta version of an automated platform with built-in SPC, intelligent analytics, and compliance-ready design.

Methods

1. Samples: Agilent MAM Sample Mix
2. Instruments: Agilent 6560
3. Platform Features:
 - Built-in SPC for LC and MS.
 - Simultaneous monitoring of multiple CQAs (e.g., glycosylation, deamidation, oxidation).
 - Real-time anomaly detection with root cause insights.
 - Visualization of historical performance data and automated reporting.
4. Comparative Analysis: Compared performance, precision, and efficiency with manual workflows.

Results and Discussion



Sample	Targets	Additional Parameters
05222025_AgilentMix_Fg100_Vcp4k_GT250_SG300_rep2	Fail	Pass
05222025_AgilentMix_Fg100_Vcp4k_GT250_SG300_rep3	Fail	Pass
05222025_AgilentMix_Fg100_Vcp4k_GT250_SG300_rep1	Fail	Pass

Product	Lot	Method	Time
Northeastern BATL SPC Fg100_Vcp4k_GT250_SG300		05222025_AgilentMix_Fg100_Vcp4k_GT250_SG300.m	01 Jan 2021 12:00:00

Results and Discussion

Benefits and Observations:

- SPC reduces downtime and variability by ensuring optimal instrument performance.
- Assessing system readiness is strengthened by a diverse set of peptides, inclusive of representative CQAs
- CQA Visualization: Enables monitoring of attribute trends for better process control.
- Dynamic results viewing to quickly assess Pass/Fail criteria
- Anomaly Detection: Flags abnormal peaks and results to streamline troubleshooting and root cause analysis, expediting investigations.

Implications:

This platform will support robust, real-time, and automated MAM implementation, optimizing product characterization and QC efficiency.

Major Findings

- Automated SPC enhances system reliability in MAM workflows.
- Compliance-ready infrastructure supports seamless QA/QC processes.
- Fragmentor voltage gradient (70V vs 100V) was used on Agilent 6560 system in lieu of skimmer voltage optimization due to instrument specifications.
- Fragmentor voltage gradient (70V vs 100V) was used on Agilent 6560 system in lieu of skimmer voltage optimization due to instrument specifications.
- Real-time data access and visualization improve decision-making.

References

1. Millán-Martín et al. (2023) *Crit Rev Anal Chem* 54:3234-3251
2. Jakes et al. (2021) *J Am Soc Mass Spectrom* 32:1998-2012
3. Yang et al. (2023) *mAbs* 15:2197668