

Developing the Research to Routine Workflows with FAIMS: Automating Large-scale SRM Method Creation for Routine HeLa Peptide Screening

Paul Humphrey,¹ Kerry Hassell², Mary Blackburn³, Michael Belford³, Michael Volny³, Scott Peterman³, Romain Huguet³; ¹Boundary Way, Hemel Hempstead, UK; ²265 Davidson Ave, Suite 101, Somerset, NJ 07783; ³355 River Oaks Parkway, San Jose, CA 951122

ABSTRACT

Purpose: To demonstrate a discovery to quantitation workflow for a peptide panel using FAIMS.
Methods: Used high resolution data to pick peptides for proteins and converted them to an SRM table to be analyzed on a triple quadrupole mass spectrometer.
Results: Improvement of 95% of the transitions monitored was facilitated by the new FAIMS source, which leads to lower limits of detection.

INTRODUCTION

Highly multiplexed protein panels are developed to enable routine sample screening while maintaining high throughput. The challenge of creating an analytically robust SRM method is the determination of which peptides to select per protein and the creation of the resulting SRM table for confident data acquisition. Each protein added to the target list increases the total SRM count by 9, quickly causing acquisition challenges on triple quadrupole mass spectrometers as most proteotypic peptides cluster into small hydrophobicity groups. To increase the selectivity space, we have incorporated a novel source, field asymmetric waveform ion mobility spectrometry (FAIMS) interface, for both profiling and screening to increase the selectivity metrics for an SRM method monitoring over 300 HeLa proteins in 60 minutes.

MATERIALS AND METHODS

Sample Preparation

A stock solution of Thermo Scientific™ Pierce™ HeLa Protein Digest was used for all experiments, with the injection of 200 ng of HeLa. Pierce Retention Time Calibration (PRTC) mixture was spiked in at 5 fmol/uL.

LC/MS

HeLa proteome profiling was performed using a Thermo Scientific™ Easy-nLC™ 1200, Thermo Scientific™ Orbitrap Tribrid mass spectrometer with a Thermo Scientific™ FAIMS™ Pro interface. A HeLa digest was injected and analyzed using a single compensation voltage (CV) setting by standard DDA methods and repeated for eight different CV settings. Each RAW file was processed to create a data matrix of proteins and peptides, retention time, CV, and precursor and product ion distribution profiles.

Data Analysis

A routine was created to construct a scheduled SRM table for the top 300 HeLa proteins using over 2,500 SRM transitions. The SRM table was imported into a triple quadrupole mass spectrometer, the Thermo Scientific™ TSQ Altis™ mass spectrometer with the FAIMS Pro interface and evaluated for analytical performance.

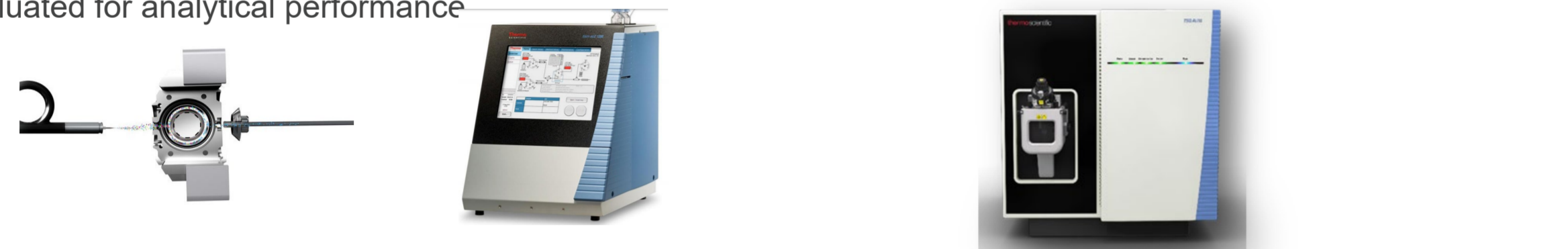


Figure 1. From left to right, the new FAIMS Pro interface, Easy-nLC 1200 system, and TSQ Altis triple quadrupole mass spectrometer.

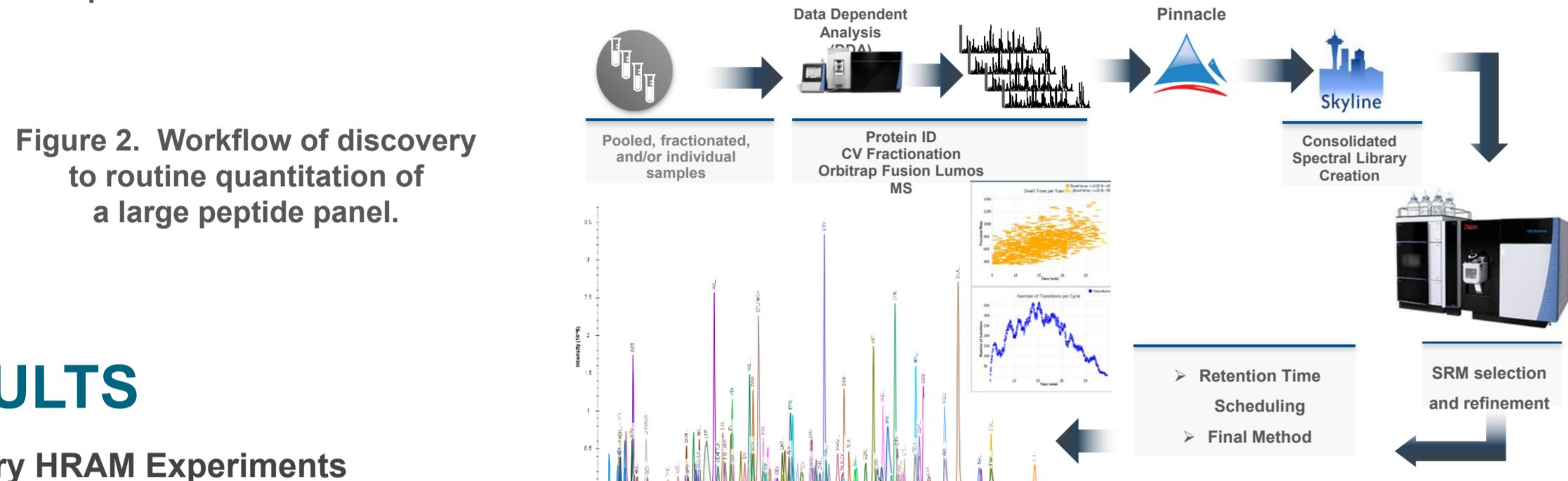


Figure 2. Workflow of discovery to routine quantification of a large peptide panel.

RESULTS

Discovery HRAM Experiments

The discovery method was used to fully characterize the HeLa digest. Replicate sample injections using single CV settings significantly increase the protein coverage from 310 proteins *without* FAIMS to over 500 proteins *with* FAIMS. Example of the base peak is shown in Figure 3, demonstrating the improvement using FAIMS.

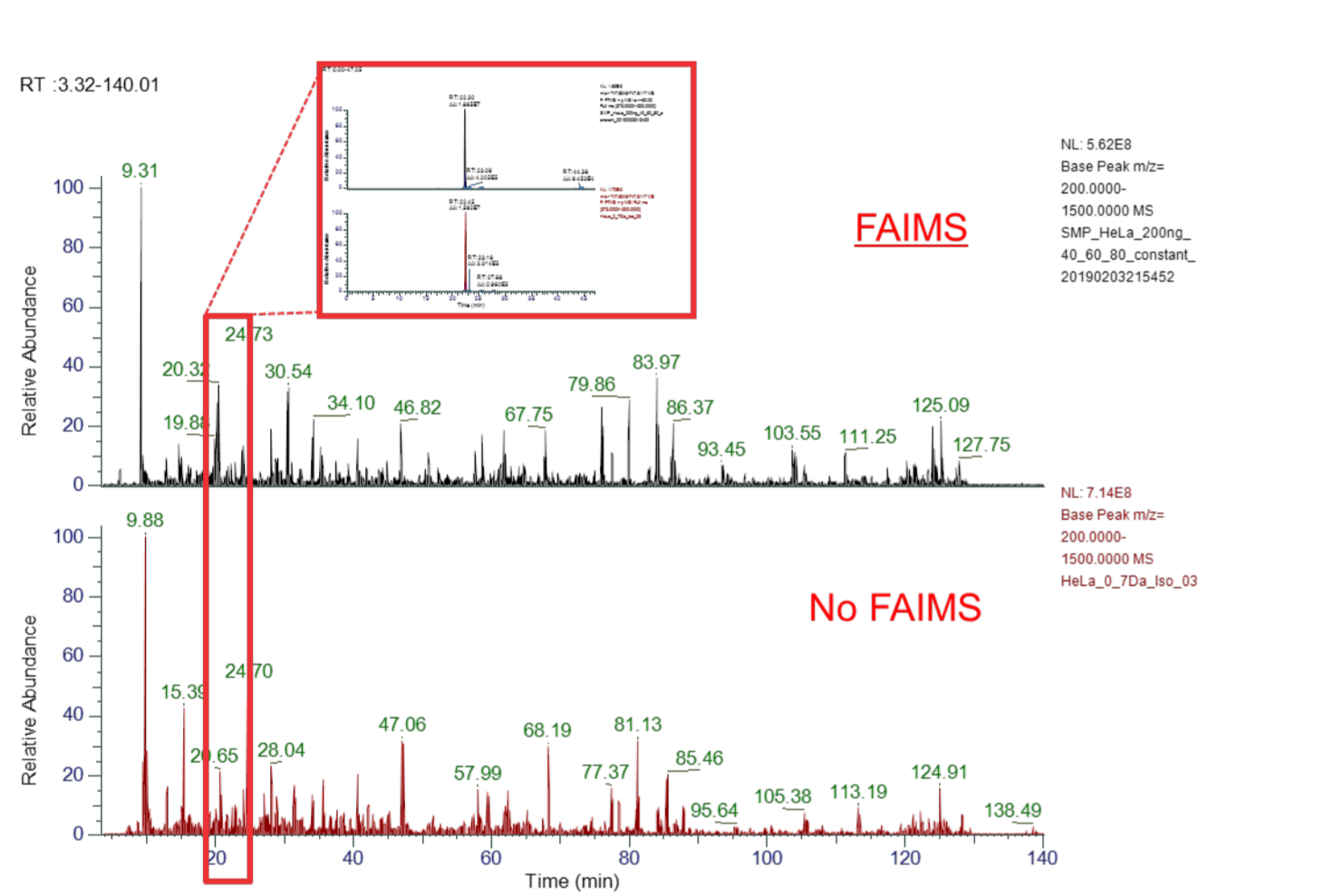


Figure 3. HRAM spectra of the overall base peak chromatogram with and without FAIMS. The inset demonstrates a low level peptide gaining intensity with FAIMS

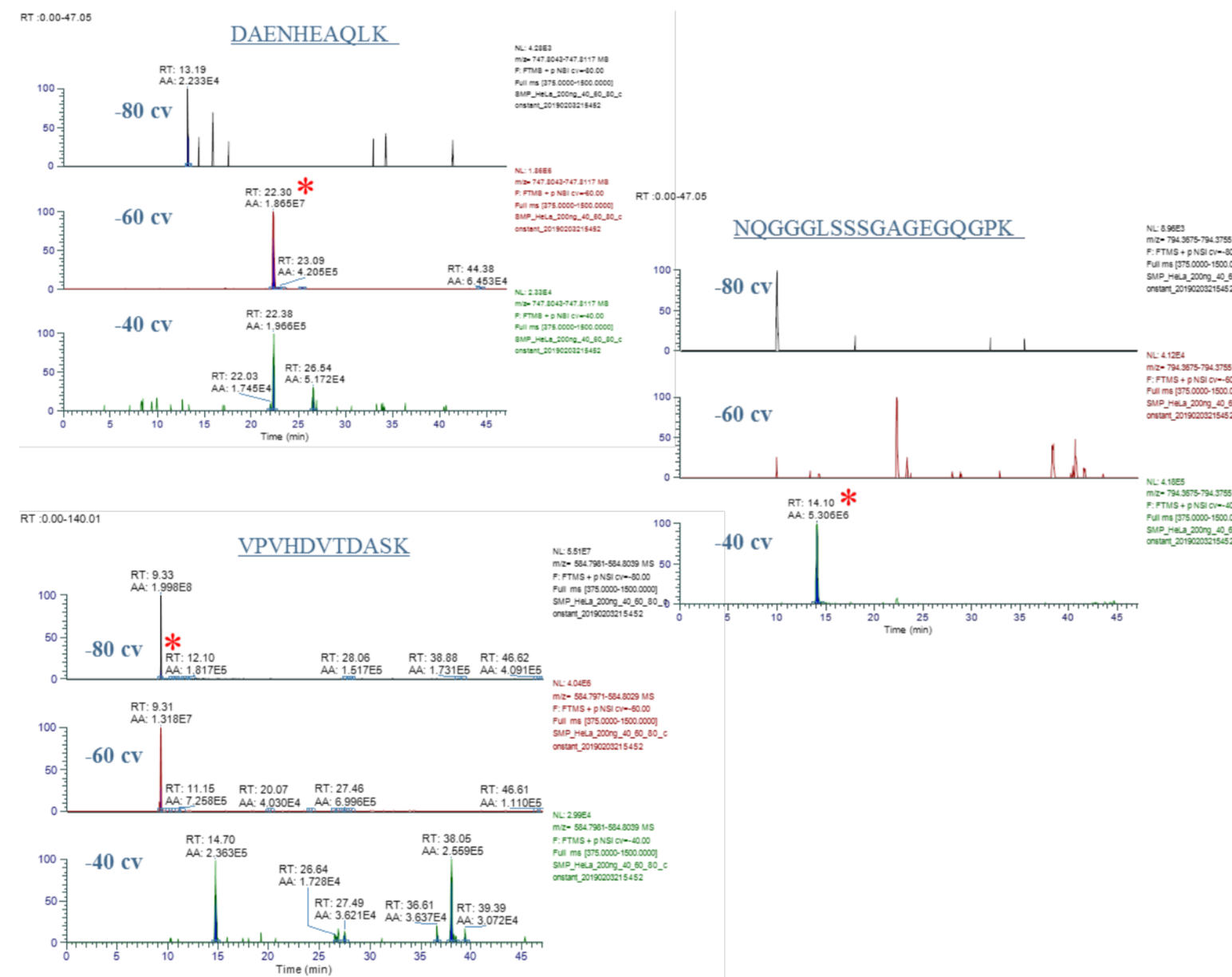


Figure 4. Showing the importance of selecting the right compensation voltage for three different peptides. (*) representing the optimal CV.

Optimization of compensation voltages is demonstrated in Figure 4 for three different peptides, showing the importance of using the correct CV for individual peptides.

Figure 5 is an example of the improvement of peptide fragmentation when using FAIMS.

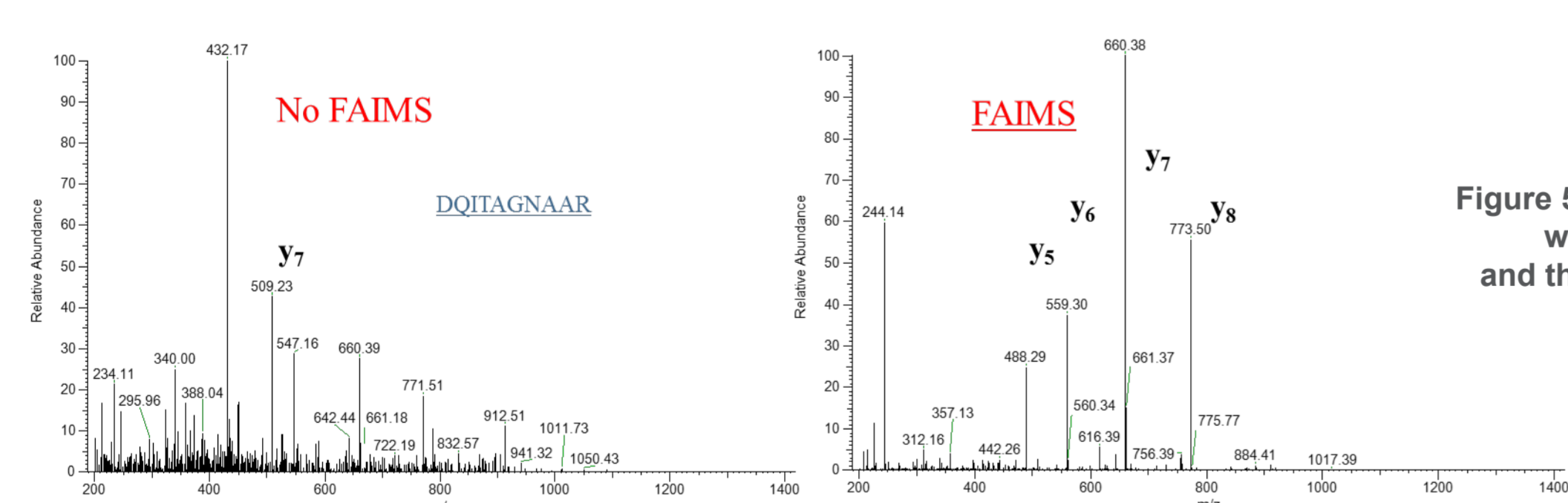


Figure 5. Spectra of peptide DQITAGNAAR without FAIMS and with FAIMS, and the respective -y ion fragmentation series of the peptides.

Peptide/Protein Selection - Pinnacle Software

Pinnacle software (Optys Tech Corporation) offers capabilities of searching routing (either spectral or sequence matching), new library creation, facilitating user-defined protein selections, establishment of peptide selection rules, and building the PRM/SRM assay. Example of the peptide selection is shown in Figure 6.

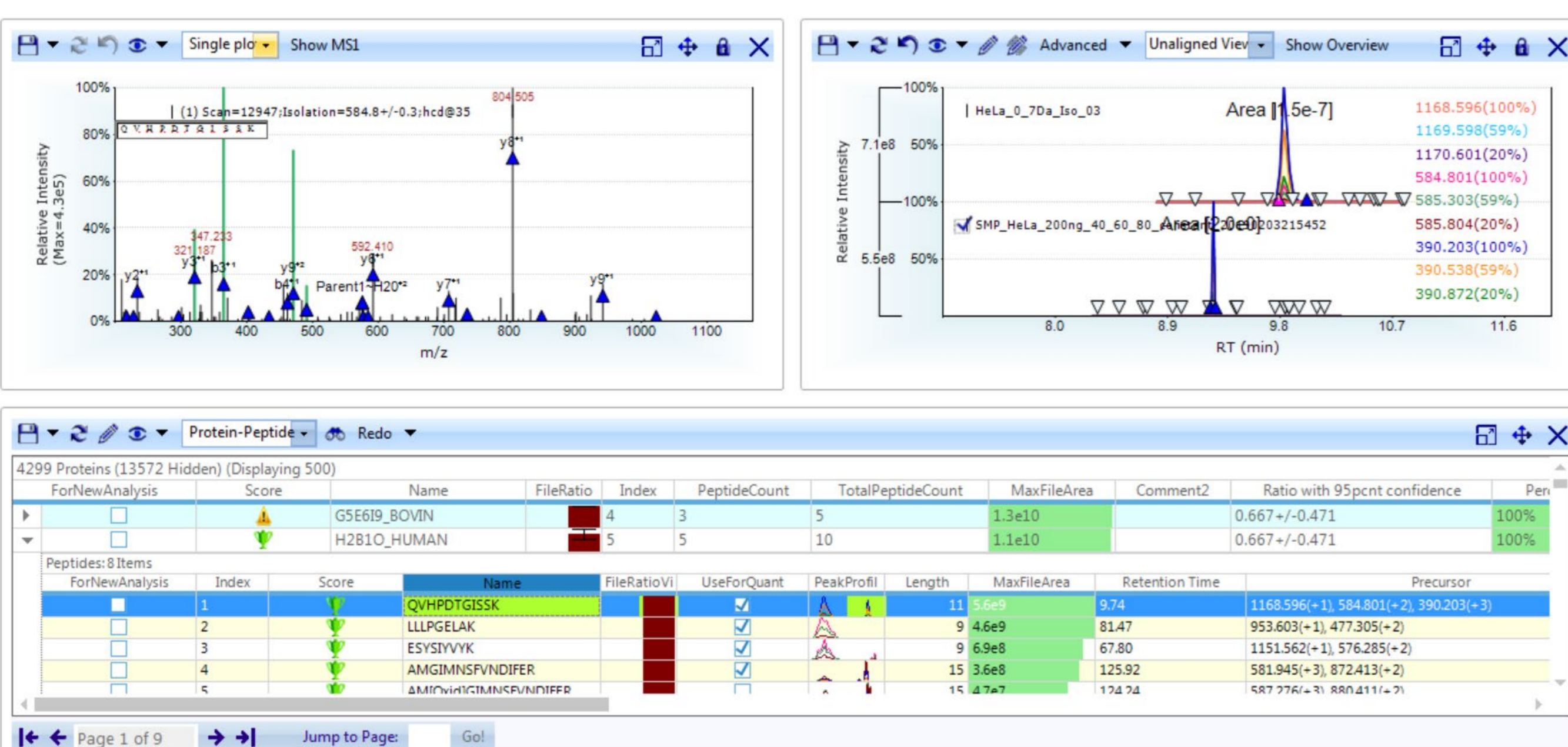


Figure 6. Screenshot of workflow of the selection of proteins/peptides from Pinnacle.

In addition, the resulting data are used to create a four-dimensional library that consists of the protein and corresponding peptides, and for each peptide, the measured retention time, CV setting, and optimal precursor m/z value and product ion distribution.

Targeted Experiments – Triple Quadrupole Mass Spectrometry

The addition of FAIMS enhances the selectivity and sensitivity of peptides, thus increasing the number of available peptides per targeted protein and resulting in more options to be considered in creating the scheduled SRM table.

The set of experiments was performed on a high-end triple quadrupole mass spectrometer capable of acquiring robust data with less than 5 msec dwell times per SRM transition and variable dwell time settings per SRM transition. The SRM table is presented in Figure 7. For each peptide, one to three transitions were selected to monitor; thus resulting in many transitions for monitoring. In order to determine if there was enough dwell time, a visualization tool in the method editor software was used to schedule windows, and is shown in Figure 8.

Compound	Retention Time (min)	RT Window (min)	Precursor (m/z)	Product (m/z)	Collision Energy (V)	Min Dwell Time (ms)	FAIMS CV (V)	
438	LAQMQDQK(-2)	24.3	496.229	876.367	19.2	10.000	-80	
439	MEDEMR(-2)	24.3	496.726	876.404	19.2	10.000	-80	
400	EAQR(-17.0214)	24.37	496.726	577.276	19.2	10.000	-80	
401	EAQR(-17.0214)	24.37	496.726	664.308	19.2	10.000	-80	
402	DHSGVAQVR(-2)	25.06	1.5	484.749	567.252	18.8	10.000	-50
404	DHSGVAQVR(-2)	25.06	1.5	484.749	794.379	18.8	10.000	-50
405	TC(-17.021460VA)	25.27	1.5	732.298	1183.442	27.2	10.000	-50
406	IAEMETQ(-2)	25.44	1.5	475.236	565.262	18.5	10.000	-50
407	IAEMETQ(-2)	25.44	1.5	475.236	636.302	18.5	10.000	-50
408	PMHEHQGQK	26.08	1.5	493.251	460.733	16.1	10.000	-80
409	PMHEHQGQK	26.08	1.5	493.251	679.316	16.1	10.000	-80
470	EEEGEEDEN(-2)	28.59	1.5	732.322	1284.449	27.2	10.000	-55
471	EEEGEEDEN(-2)	28.59	1.5	732.322	1285.553	27.2	10.000	-55
472	HLHAQSTLDAN(-)	29	1.5	661.346	970.495	24.8	10.000	-50
473	VDC(-17.021464H)	29.25	1.5	699.292	1137.43	26.1	10.000	-60
474	VDC(-17.021464H)	29.25	1.5	699.292	1183.483	26.1	10.000	-60
475	HPQQEELK(-2)	29.31	1.5	536.275	604.33	20.5	10.000	-80
476	HPQQEELK(-2)	29.31	1.5	536.275	725.321	20.5	10.000	-80
477	ARQGGTSPHLGQK	29.44	1.5	561.285	793.42	18.1	10.000	-80
478	ARQGGTSPHLGQK	29.44	1.5	561.285	1067.538	18.1	10.000	-80

Figure 7. SRM table of peptides targeted in HeLa analysis. Optimized CVs are represented in the far right column.

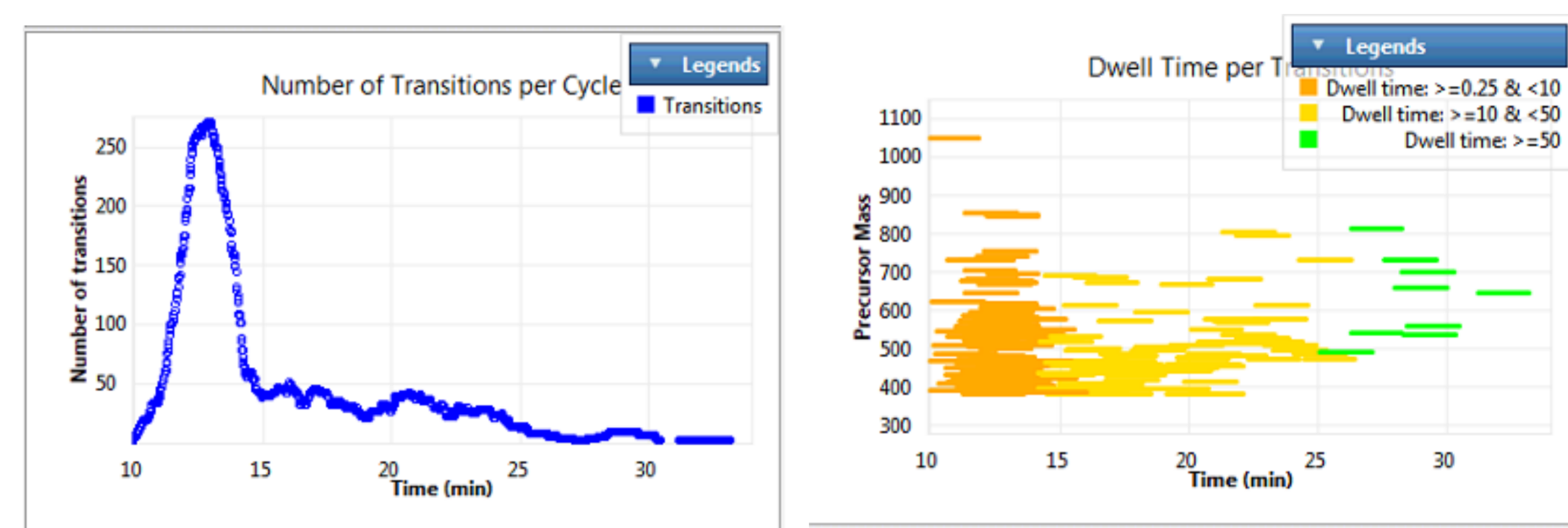


Figure 8. Visualization tool with the method editor showing the number of transitions per time, as well as the dwell times of different precursor masses over the chromatographic timescale.

For the different protein groups targeted, the optimal peptides were selected based on relative response in the discovery method, but more importantly on the retention time and CV setting as the two values were used to create the final SRM table. Peptides were grouped into overlapping retention time and CV bins to maximize duty cycle while maintaining analytical performance. Figure 9 depicts the variation of transmitted ion abundance with CV value for two peptides.

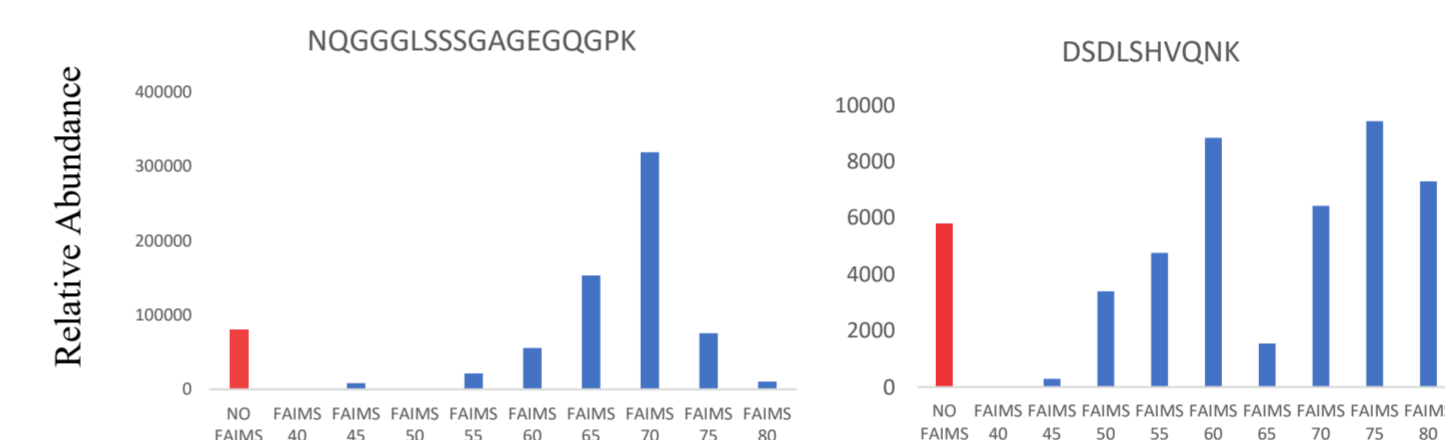


Figure 9. Comparison of peak area of no FAIMS (red) to different compensation voltages highlighted in blue.

PRTC, a well-known heavy labeled peptide standard was used to show overall intensity improvement using FAIMS, and the results can be seen in Figure 10. The standard was then made into a dilution series to study if linearity or lower limit of detection can be obtained using FAIMS. Figure 11 illustrates that the linearity is not affected by FAIMS, and Figure 12 demonstrates the ability to reach a lower LLOQ because of diminishing interferences in the quadrupole isolation window.

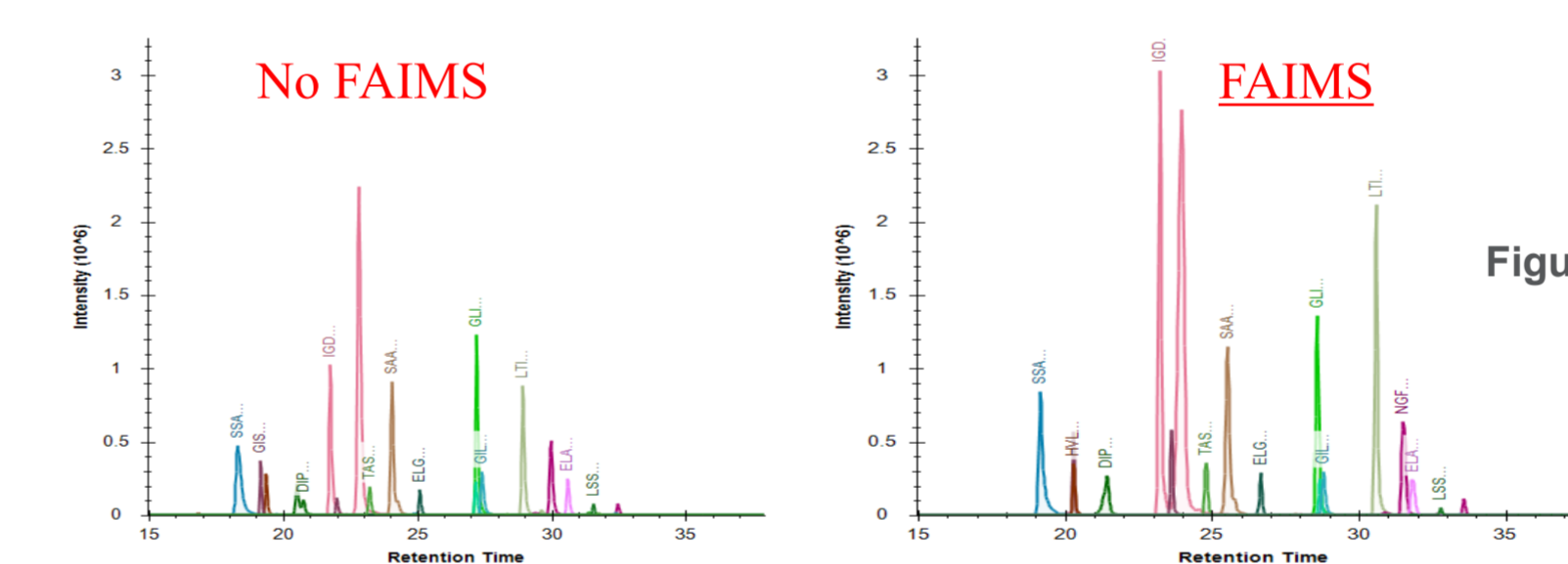


Figure 10. PRTC total ion current of peak intensity with and without FAIMS.

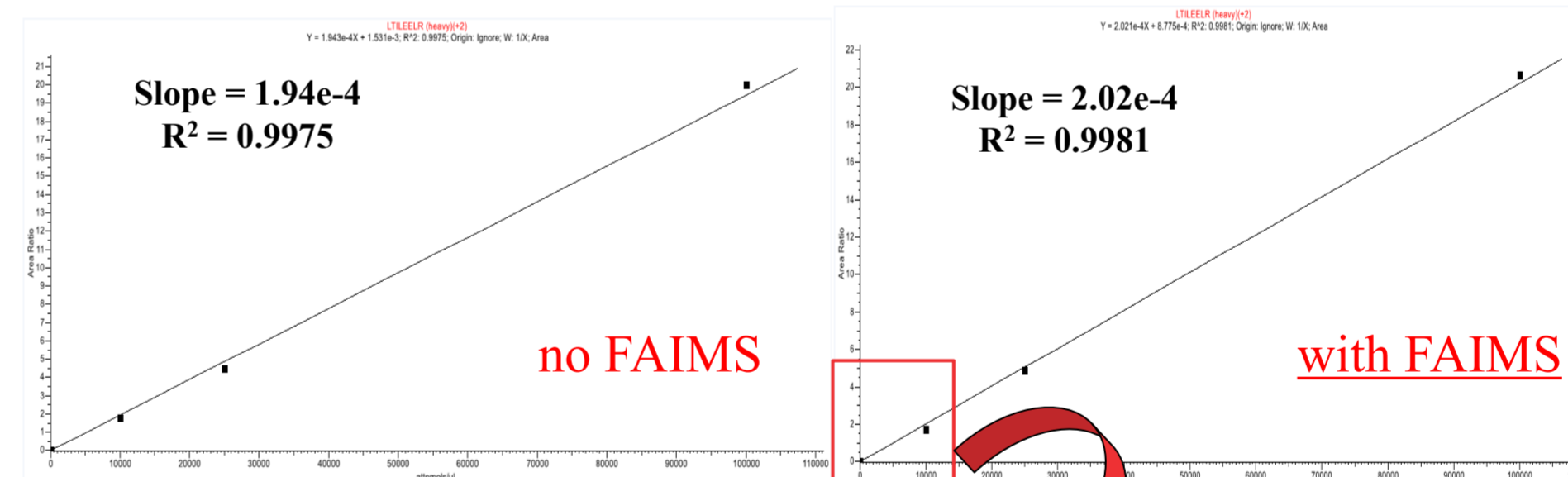


Figure 11. Peptide LTILEELR (heavy labeled) spiked into HeLa in concentration range 1 attomol/uL – 100 fmol/uL linearity remains consistent with FAIMS across linear range.

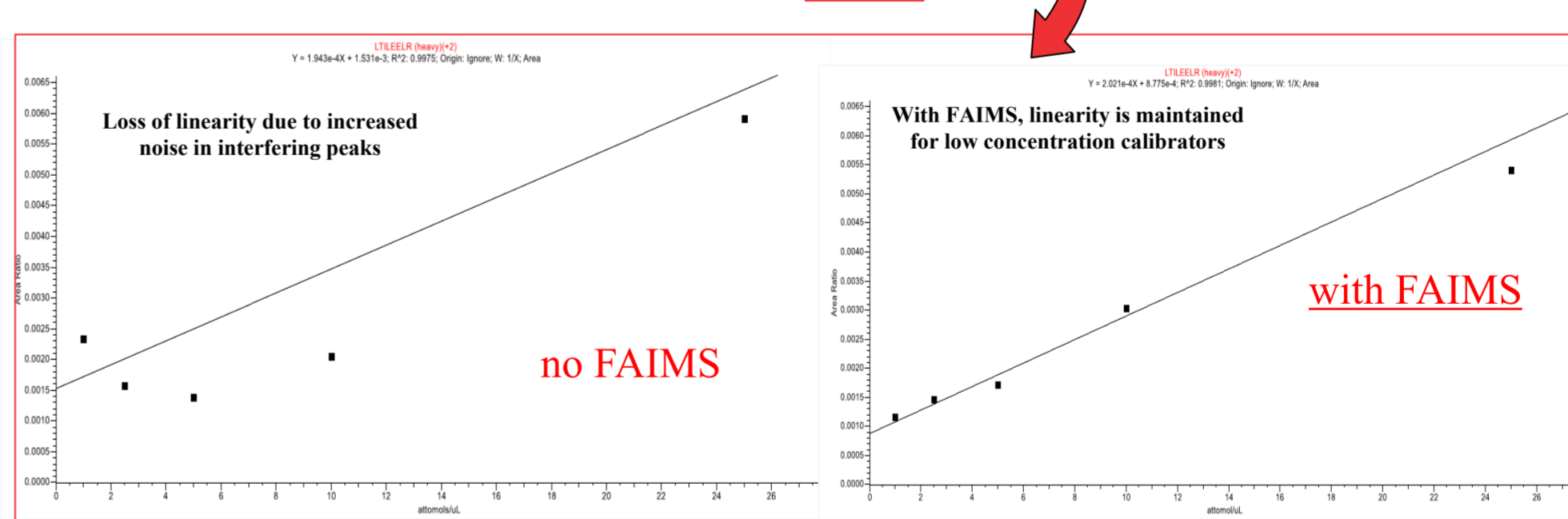


Figure 12. Peptide LTILEELR (heavy labeled) spiked into HeLa in concentration range 1 attomol/uL – 1 fmol/uL at low concentration interferences are present and with FAIMS the linearity is improved at lower concentrations.

In the 150 proteins monitored, there was a 95% improvement of peptide signal, with varying results of improvement.

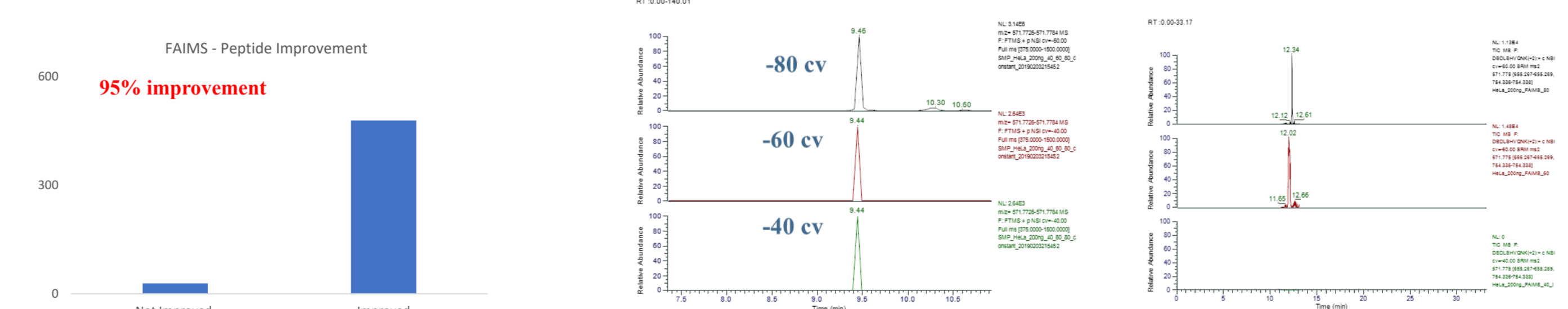


Figure 13. Consistency of source and instrumentation. The figure to the left is HRAM data at different compensation voltage, whereas the right figure is compensation voltages on a triple quadrupole. The same CV is ideal for this peptide on both instruments

CONCLUSIONS

- Field asymmetric waveform ion mobility spectrometry (FAIMS) can be used in a discovery environment to determine a peptide screening panel. These peptides can be put into a targeted panel to be monitored by a triple quadrupole mass spectrometer. The use of the FAIMS Pro interface increases signal-to-noise of these peptides.
- Different compensation voltages can be run early in the workflow process to determine the optimal voltage this is then translated to the targeted panel without additional optimization.
- Due to the improvement of signal-to-noise, linearity of calibration curves is improved and lower LLOQs can be obtained.
- Further analysis will be done to inquire on looking at the number of transitions per peptide and determining what protein expression levels can be monitored

For Research Use Only. Not for use in diagnostic procedures.

TRADEMARKS/LICENSING

© 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.