

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

Kerry Hassell¹, Mary Blackburn², Michael Belford², Michael Volny², Scott Peterman², Romain Huguet²; ¹ 265 Davidson Ave, Suite 101, Somerset, NJ 07783, ² 355 River Oaks Parkway, San Jose, CA 95112

ABSTRACT

Purpose: To demonstrate a discovery to quantitation work flow for 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 seen using the new FAIMS source which lead to lower limits of detection.

INTRODUCTION

Highly multiplexed protein panels are developed to enable routine sample screening while maintaining high throughput. The challenge to creating an analytically robust SRM method is determining which peptides to select per protein and creating the resulting SRM table for confident data acquisition. Each protein added to the target list increases 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, injection 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 an 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 2500 SRM transitions. The SRM table was imported into a triple quadrupole mass spectrometer, Thermo Scientific™ TSQ Altis™ with the FAIMS Pro interface and evaluated for analytical performance.

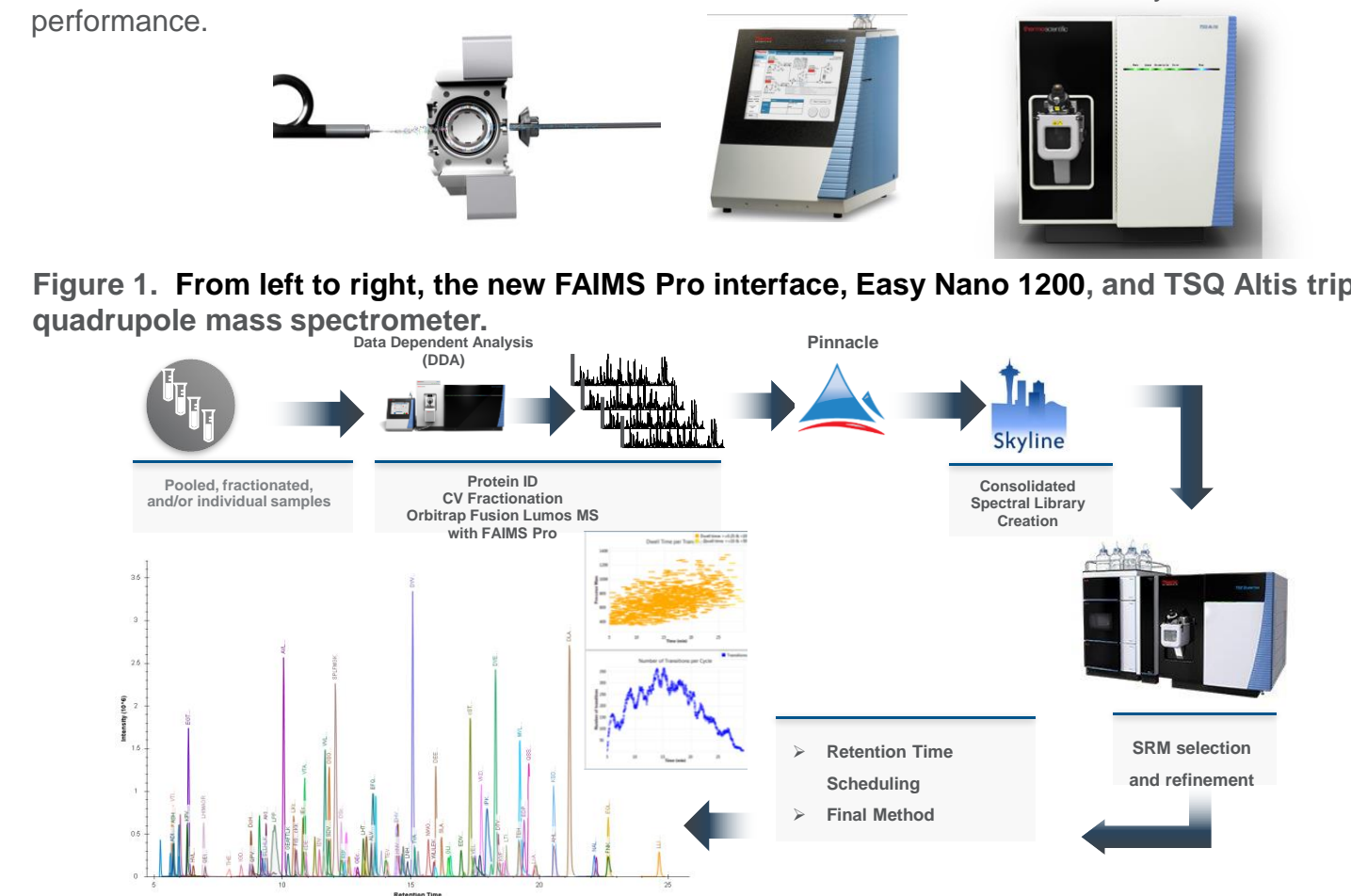


Figure 2. Workflow of discovery to routine quantitation of a large peptide panel.

RESULTS

Discovery HRAM Experiments

The discovery method was used to fully characterize the HeLa digest. Replicate sample injections using single CV settings significantly increases 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.

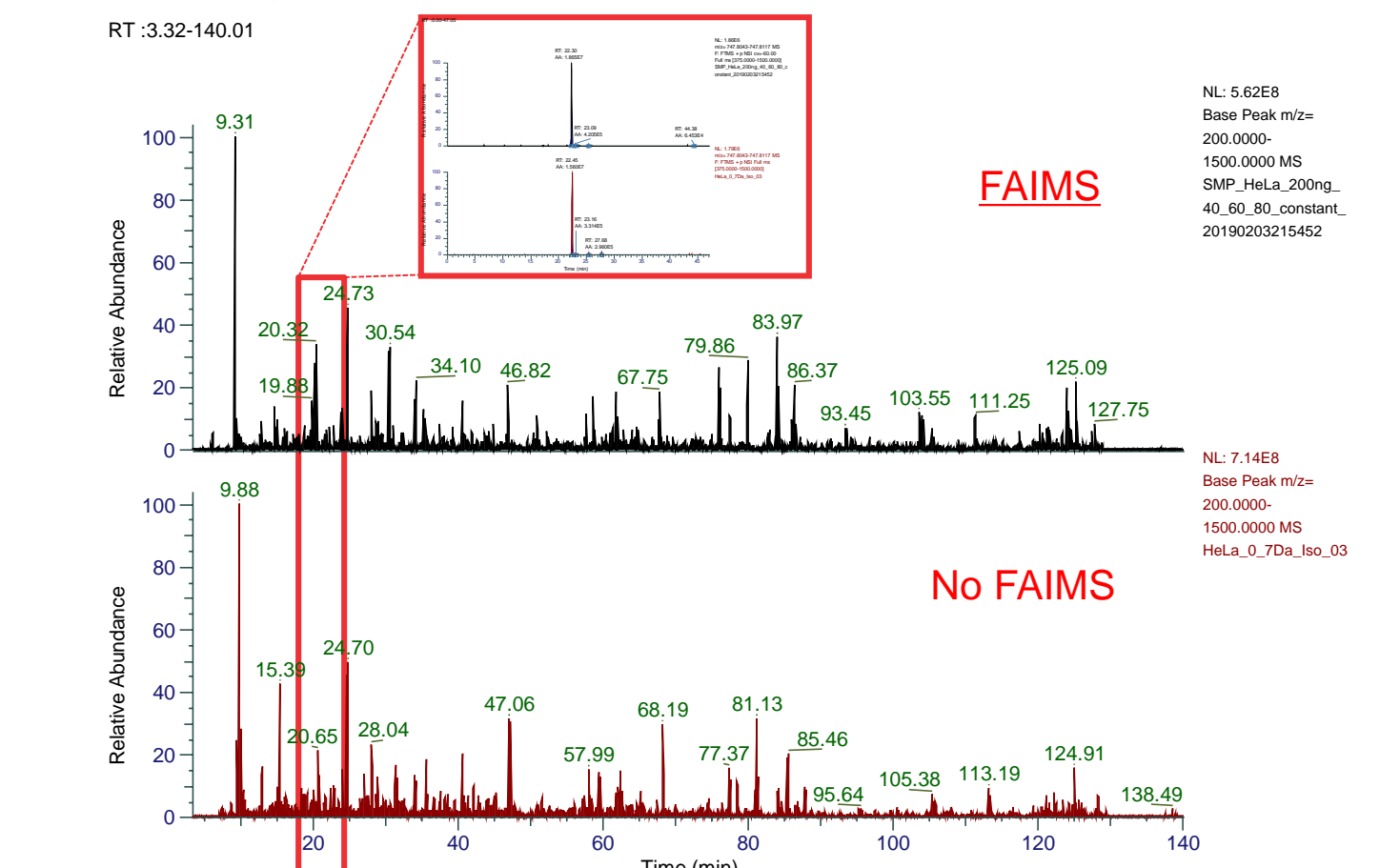
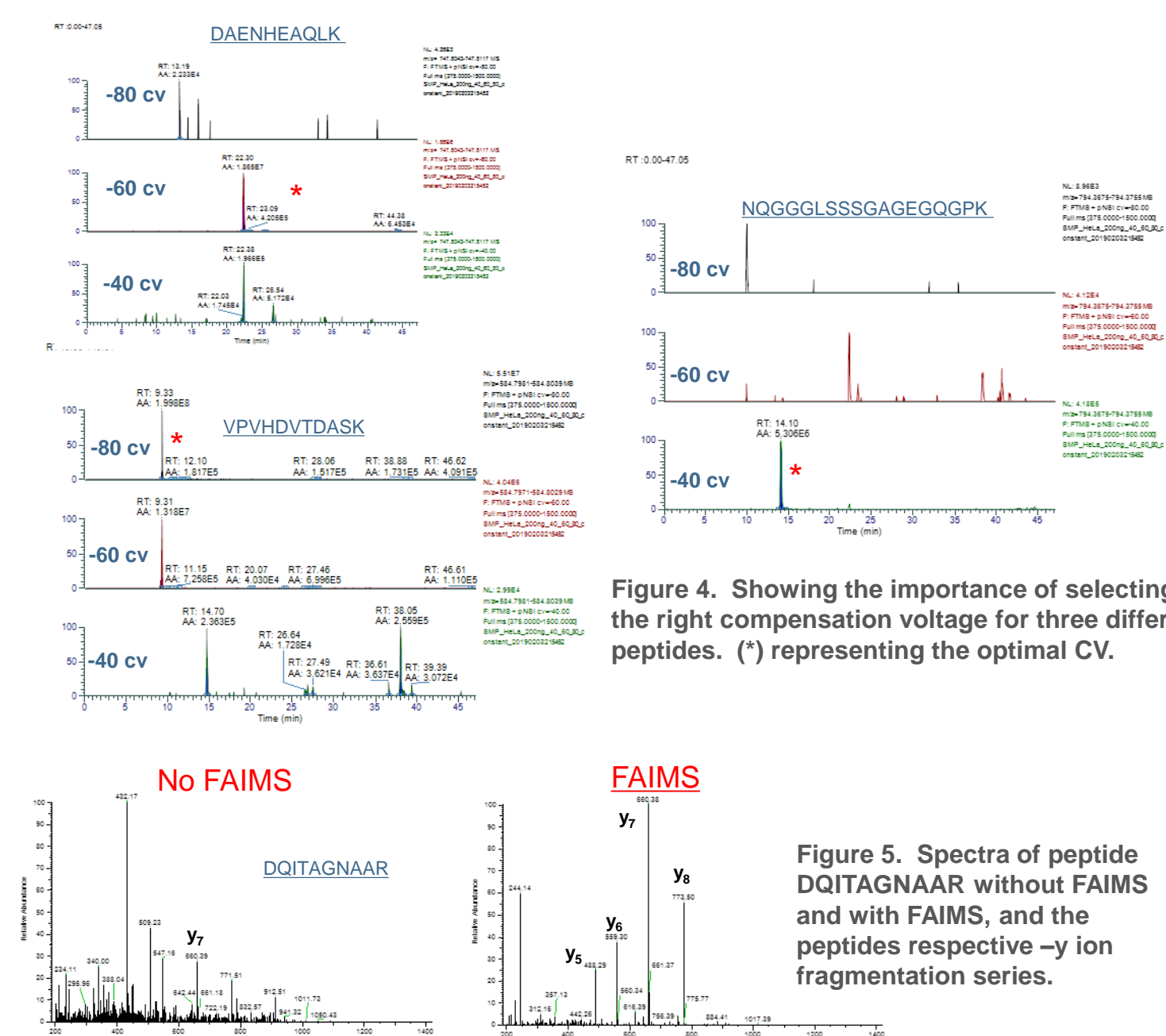


Figure 3. HRAM spectra of the overall base peak chromatogram with and without FAIMS. The inset demonstrates a low level peptide gaining intensity with FAIMS

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.



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, 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; this resulted in many transitions to be monitored. 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 (eV)	Min Dwell Time (ms)	FAIMS CV (V)
436 UNKQHCQYDQ	26.3	1.5	496.226	88.387	19.2	110.027	80
438 NEDSMRRI-2	26.3	1.5	496.726	879.484	19.2	110.027	80
440 EAQSQI-17-02241	24.37	1.5	496.726	577.276	19.2	110.027	80
441 EAQSQI-17-02241	24.37	1.5	496.726	664.308	19.2	110.027	80
442 DSPHGAVPR-2	25.06	1.5	484.749	567.252	18.8	110.027	50
443 DSPHGAVPR-2	25.06	1.5	484.749	572.351	18.8	110.027	80
444 DSPHGAVPR-2	25.06	1.5	484.749	794.579	18.8	110.027	50
445 TCY-17-0214601	29.37	1.5	732.302	120.542	27.2	110.027	80
446 IAMMTRDQ-2	25.44	1.5	473.266	305.362	18.5	110.027	80
447 IAMMTRDQ-2	25.44	1.5	473.266	636.302	18.5	110.027	80
448 IHMAINQPSQI	26.08	1.5	493.251	460.733	16.1	110.027	80
449 IHMAINQPSQI	26.08	1.5	493.251	679.304	16.1	110.027	80
470 EESAESEDSQD-2	25.59	1.5	732.322	120.449	27.2	110.027	55
471 EESAESEDSQD-2	25.59	1.5	732.322	120.553	27.2	110.027	65
472 HETHAQSLDIAK	29	1.5	661.346	970.495	24.8	110.027	50
473 VDQI-17-0214611	29.25	1.5	699.202	1137.43	26.1	110.027	40
474 VDQI-17-0214611	29.25	1.5	699.202	1183.463	26.1	110.027	40
475 NPQDQELHAI-2	29.31	1.5	536.273	604.305	20.5	110.027	40
476 NPQDQELHAI-2	29.31	1.5	536.273	723.323	20.5	110.027	40
477 HQTGDSHSLQDQI	29.44	1.5	561.285	790.642	18.1	110.027	40
478 HQTGDSHSLQDQI	29.44	1.5	561.285	1067.333	18.1	110.027	40

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

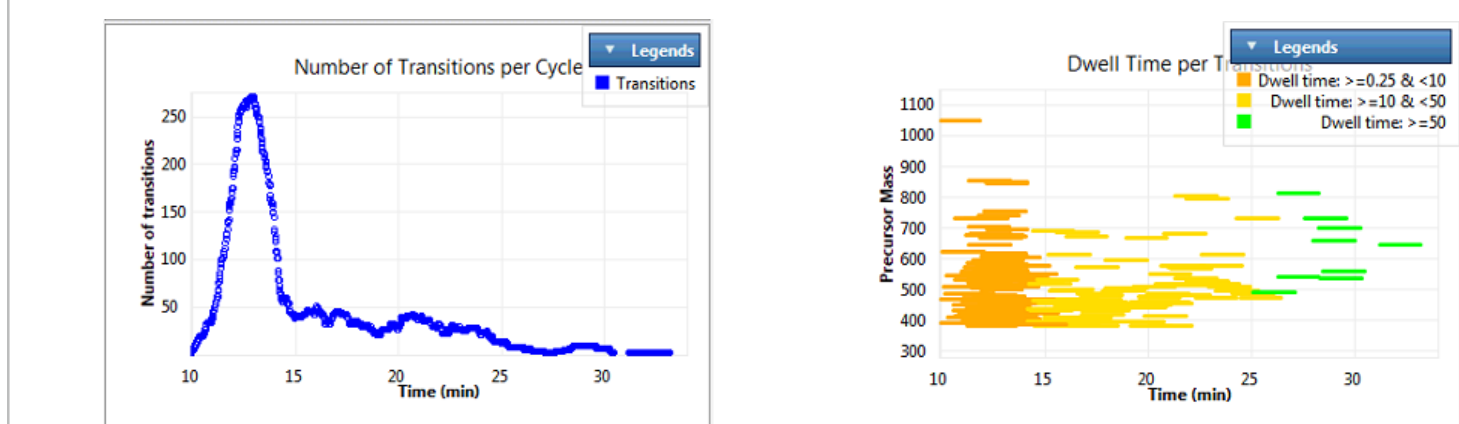


Figure 8. Visualization tool with the method editor showing the number of transitions per time as well as the dwell times of different precursor masses over the chromatographic timescale.

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. FIG. 7 depicts the variation of transmitted ion abundance with CV value for two peptides.

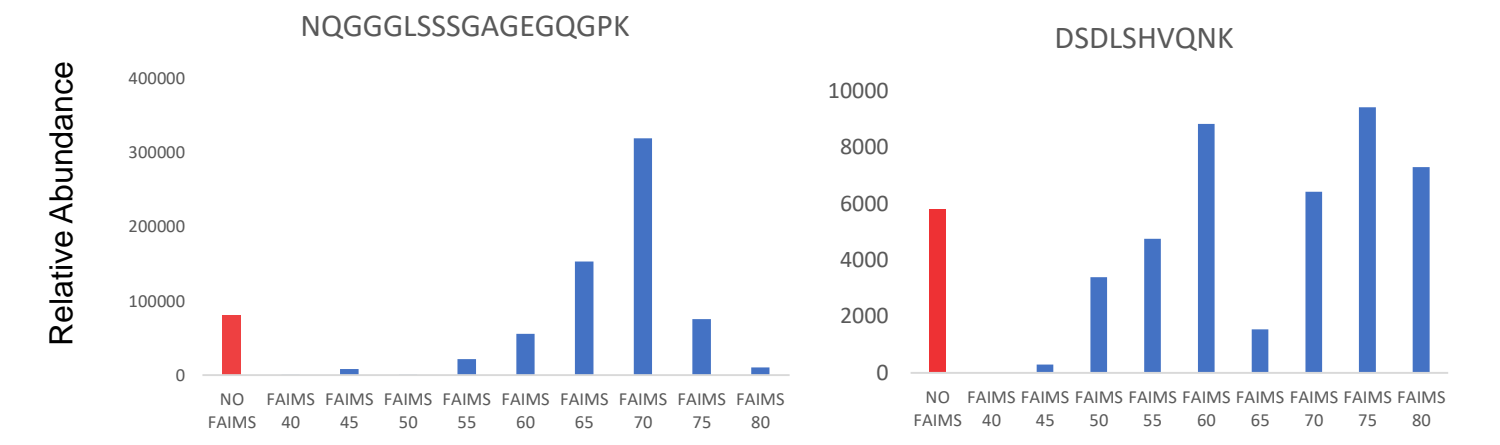


Figure 7. Comparison of peak area of no FAIMS, red, to different compensation voltages highlighted in blue.

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 8. The standard was than made into a dilution series to study if linearity or lower limit of detection can be obtained using FAIMS. Figure 8 illustrates that the linearity is not affected by FAIMS, and Figure 9 demonstrates the ability to reach a lower LLOQ because of diminishing interferences in the quadrupole's isolation window.

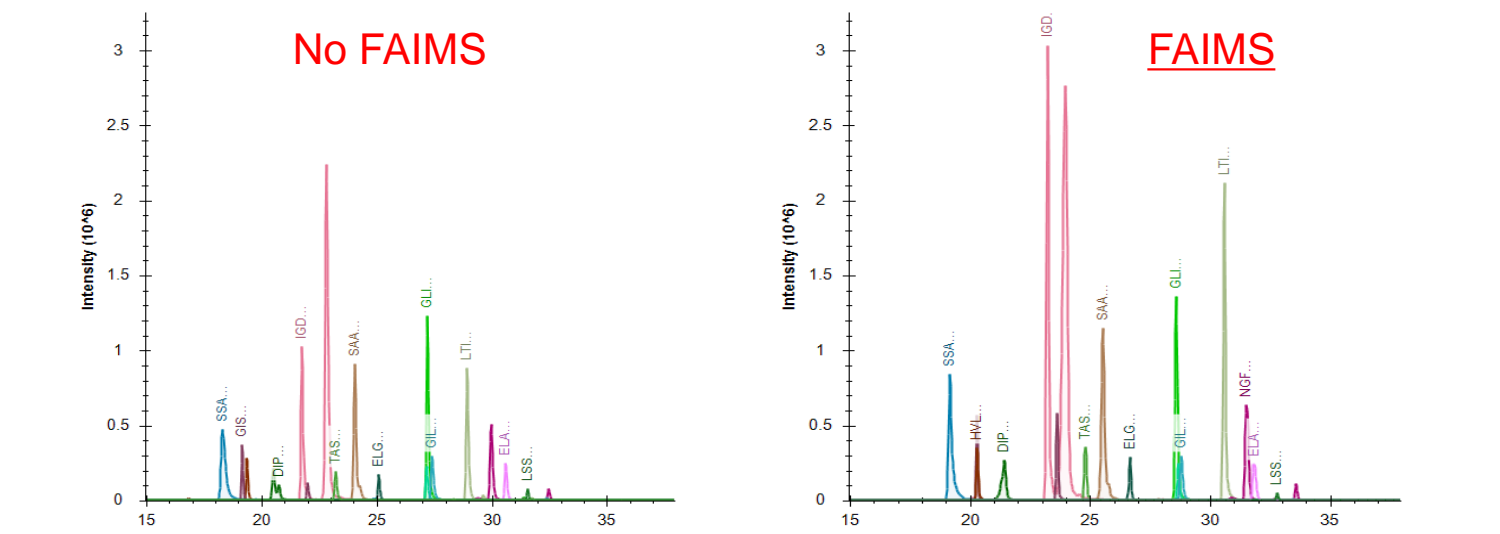


Figure 8. PRTC total ion current of peak intensity with and without FAIMS.

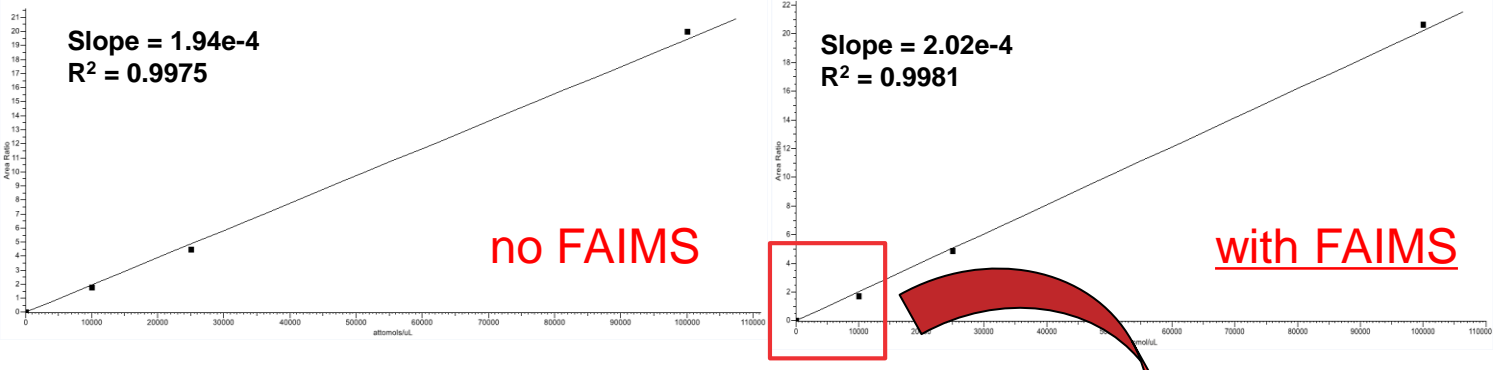


Figure 9. Peptide LTILEELR (heavy labeled) spiked into HeLa in concentration range 1 attomol/uL – 100 fmol/uL linearity remains consistent with FAIMS across linear range.

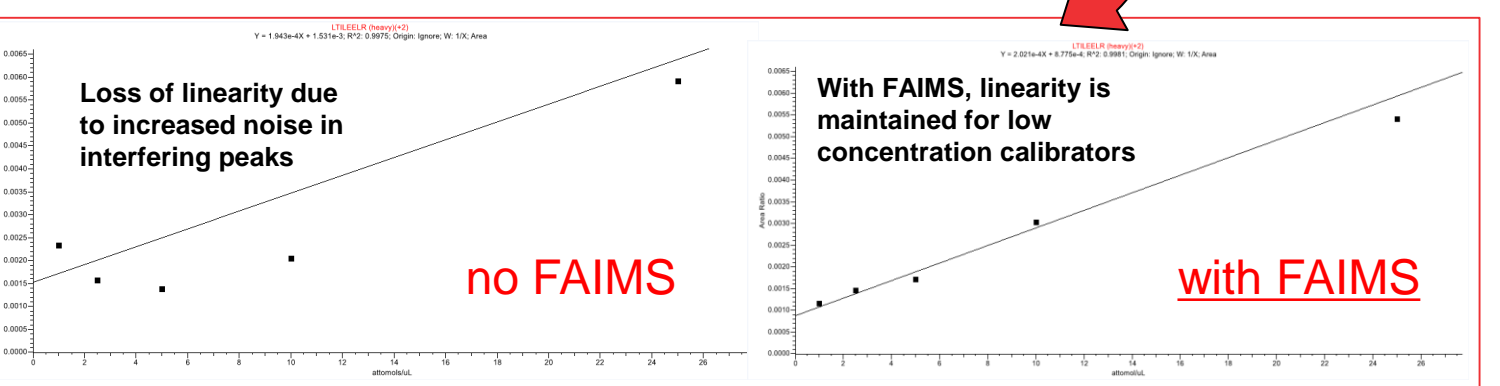


Figure 10. Peptide LTILEELR (heavy labeled) spiked into HeLa in concentration range 1 attomol/uL – 1 fmol/uL at low concentration interferences are present and with FAIMS the linearity is improved at lower concentrations.

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

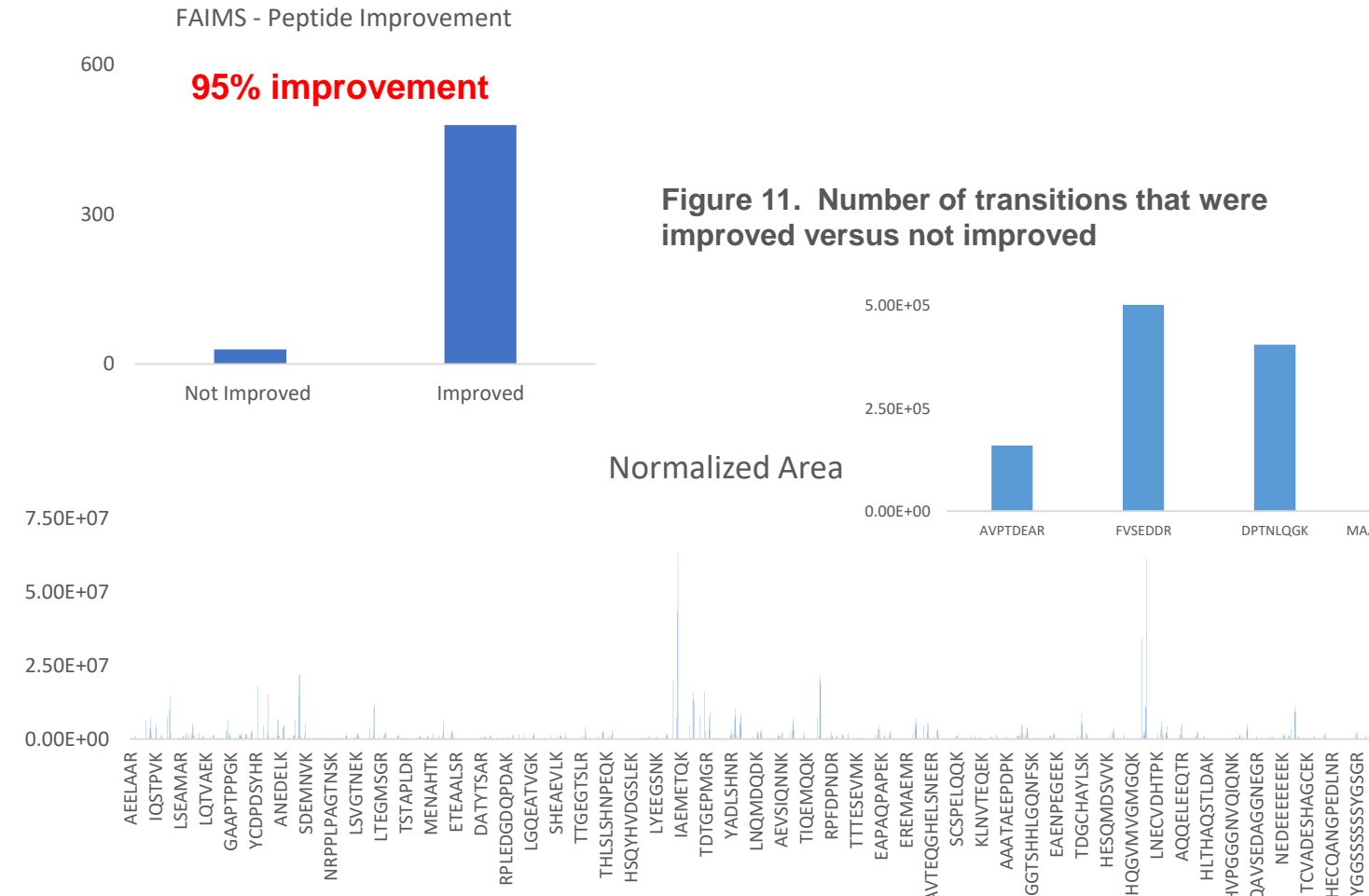


Figure 11. Number of transitions that were improved versus not improved

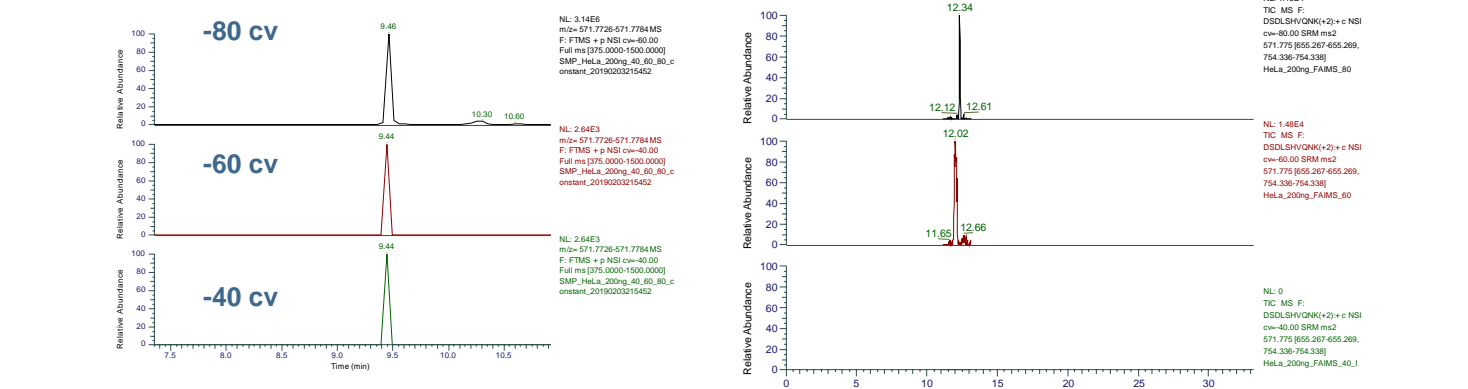


Figure 12. Consistency of source and instrumentation. The figure to the left is HRAM data at different compensation voltage, whereas the right figure is compensation voltages on a triple quadrupole. The same CV is ideal for this peptide on both instruments

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.

