

FAIMS Pro Interface Coupled to Triple Quadrupole Mass Spectrometer for Quantification of Peptides in Complex Matrices

Michael Volný, Claudia P. B. Martins, Michael W. Belford, Mary L. Blackburn, Thermo Fisher Scientific, San Jose, CA, United States

ABSTRACT

Purpose: The Thermo Scientific™ FAIMS Pro™ interface provides the selectivity and ease-of-use for improved peptide quantifications in complex matrices and for lower abundance peptides.

Methods: Thermo Scientific™ EASY-nLC™ system was used with Thermo Scientific™ TSQ Altis™ Triple Quadrupole Mass Spectrometer

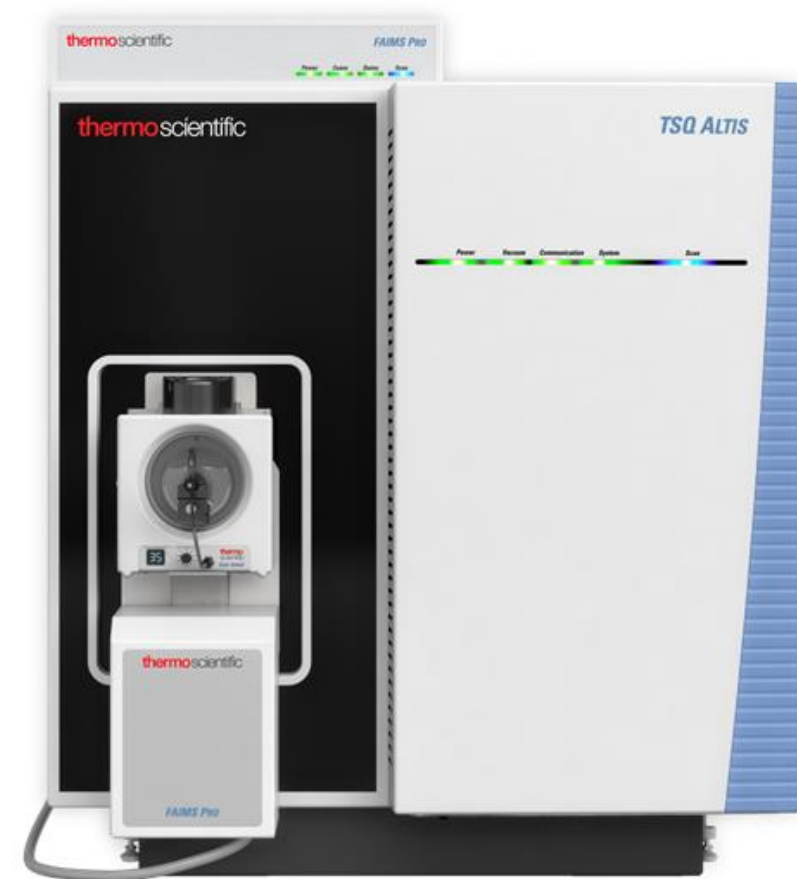
Results: Results show compatibility of FAIMS Pro interface with TSQ Altis Triple Quadrupole Mass Spectrometer and the applicability of this coupling for improvement of current peptide quantification workflows.

INTRODUCTION

The FAIMS Pro interface provides orthogonal precursor ion selectivity based on differential gas phase mobility. The Compensation Voltage (CV) setting determines which groups of ions are transmitted to the mass spectrometer for detection. A wide range of possible CV settings increases instrument performance for proteomic experiments. The increased selectivity and sensitivity enables researchers to improve quantification and to maximize efficiency and conserve sample. Many important peptides are biologically highly active and thus need to be monitored and quantified at ultra-low concentrations. The task of targeted peptide quantification is typically executed by LC-MS using triple quadrupoles in SRM/MRM mode. Such experiment can be disrupted by high background noise or interference from isobaric species. The application of differential mobility, enabled by the FAIMS Pro interface, offers an orthogonal mode of separation that enhances specificity and reduces the background interference resulting in improved LoQ levels.

MATERIALS AND METHODS

A TSQ Altis Triple Quadrupole Mass Spectrometer coupled with a Thermo Scientific™ EASY-nLC™ system was utilized for all measurements. Peptide standards and test matrices were obtained from Bachem, Thermo Fisher Scientific and Sigma-Aldrich. The FAIMS Pro interface attaches to the mass spectrometer with a flange and can be installed with Thermo Scientific NG sources.



Data Analysis

Thermo Scientific™ TraceFinder™ Software, Thermo Scientific™ FreeStyle™ Software and Skyline were used to process the results

FAIMS Pro software implementation

No Instrument configuration necessary, FAIMS Pro interface is recognized by the software when mounted and powered up

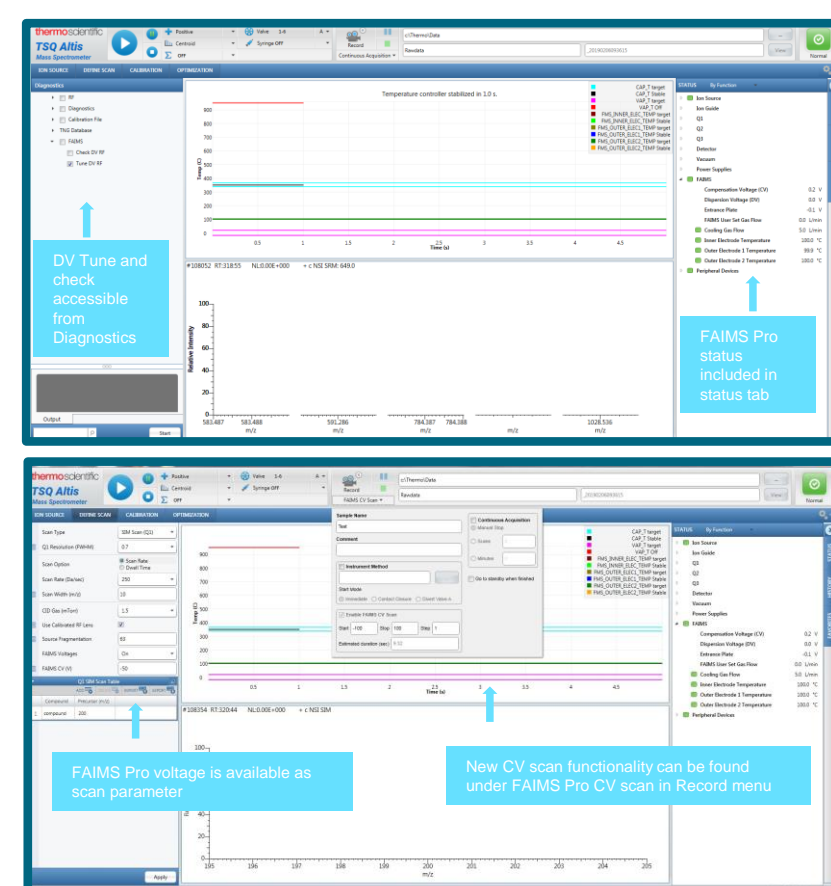
Initial DV Tune is available from Tune GUI

FAIMS parameters become part of Source and Scan tabs

CV optimization is available in Tune

Optimization files are automatically saved for further interrogation

CV value can be part of SRM or SIM table in the LC-MS method for each precursor ion



RESULTS

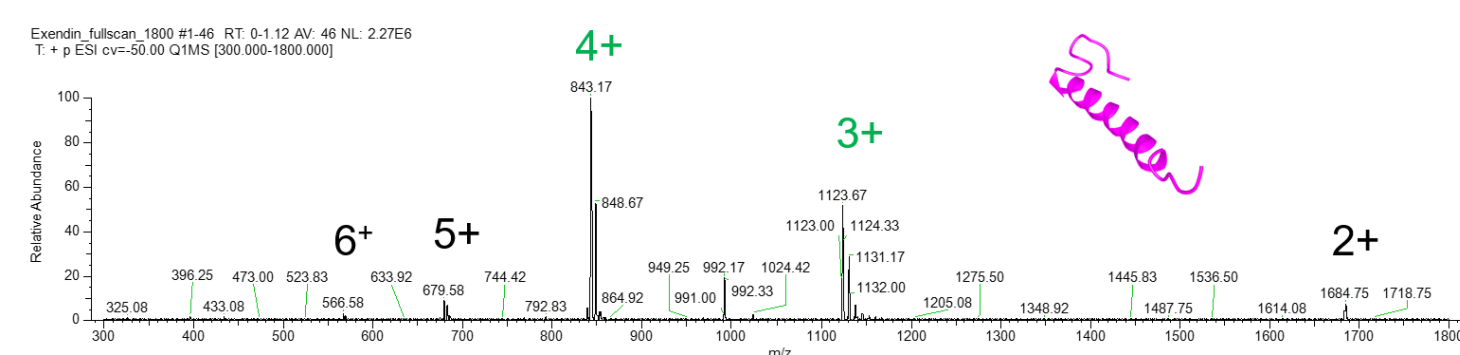


Figure 1. [9-39]Exendin (DLSKMEEAEVRLFIWELKNGPSSGAPPPS – NH₂) direct infusion ESI-MS spectrum showing the charge state distribution

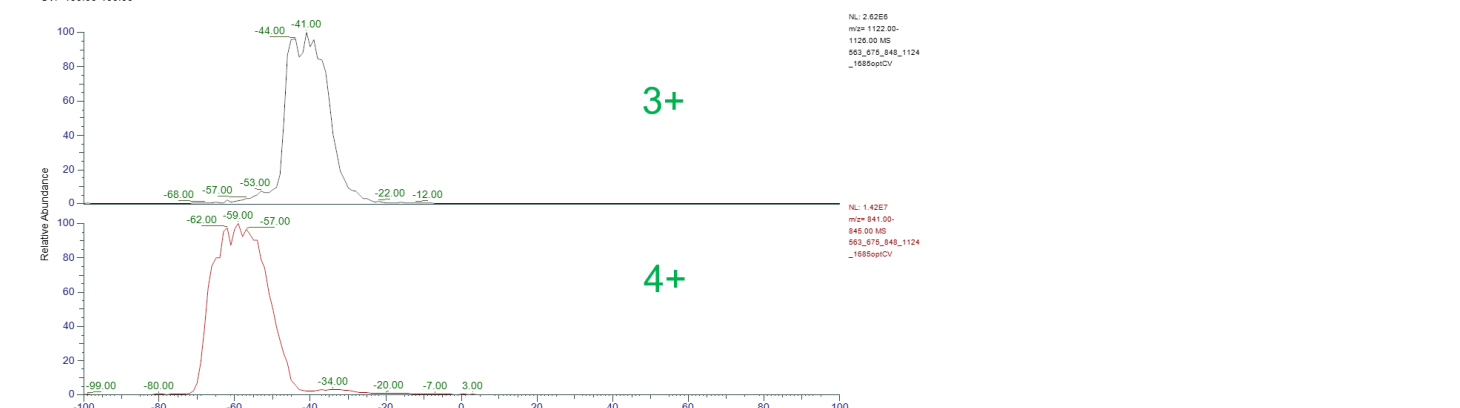


Figure 2. FAIMS Pro Compensation Voltage (CV) scan showing separation of Exendin [M+3H]⁺ and [M+4H]⁺ ions by FAIMS Pro

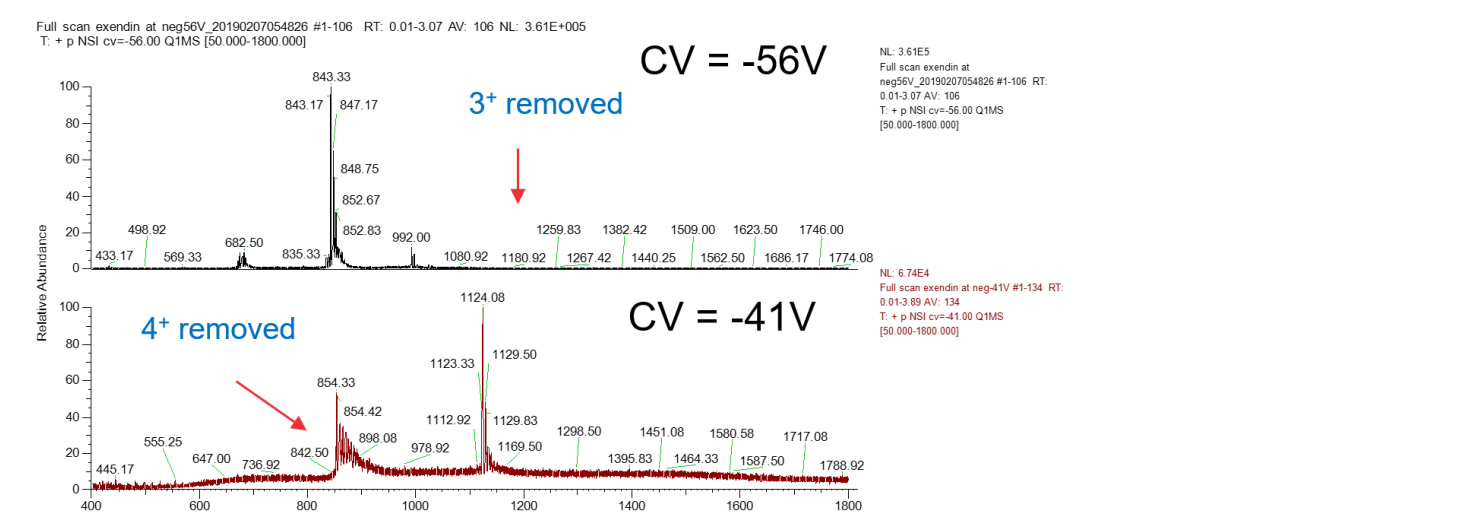


Figure 3. Setting CV to a specific value allows selective transmission of targeted charge state with removal or significant suppression of other charge states

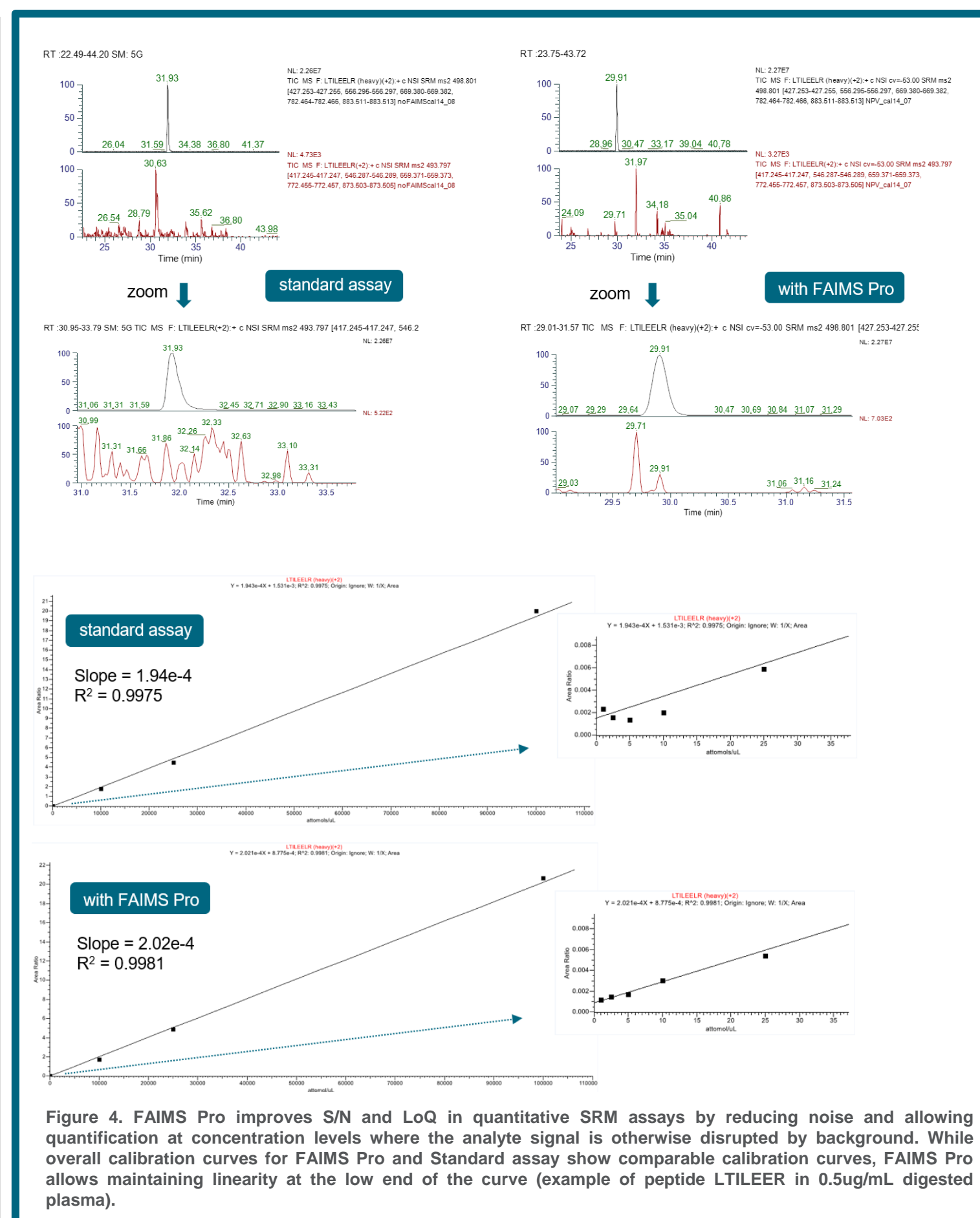


Figure 4. FAIMS Pro improves S/N and LoQ in quantitative SRM assays by reducing noise and allowing quantification at concentration levels where the analyte signal is otherwise disrupted by background. While overall calibration curves for FAIMS Pro and Standard assay show comparable calibration curves, FAIMS Pro allows maintaining linearity at the low end of the curve (example of peptide LTILEELR in 0.5ug/mL digested plasma).

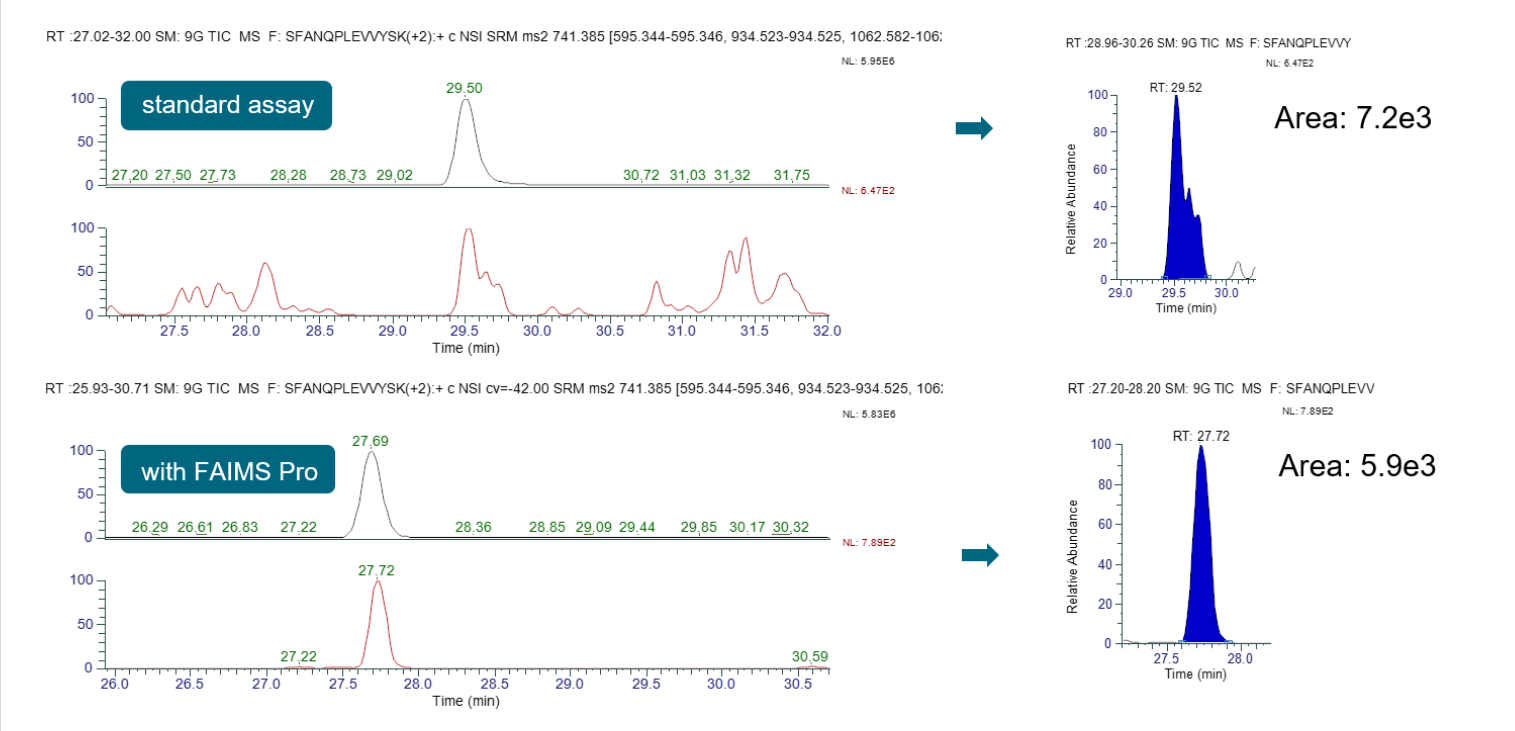


Figure 5. Example of peptide SFANQPLEVYV in digested plasma matrix shows how FAIMS Pro (bottom panel) improves peak shape by removing a co-eluting interference. As a result, automatic integration provides more accurate value without a necessity to reintegrate manually as would be required in the standard assay (top panel)

Thermo Scientific™ Pierce™ LC-MS/MS System Suitability Standard (7 x 5 Mix) with FAIMS Pro on TSQ Altis MS

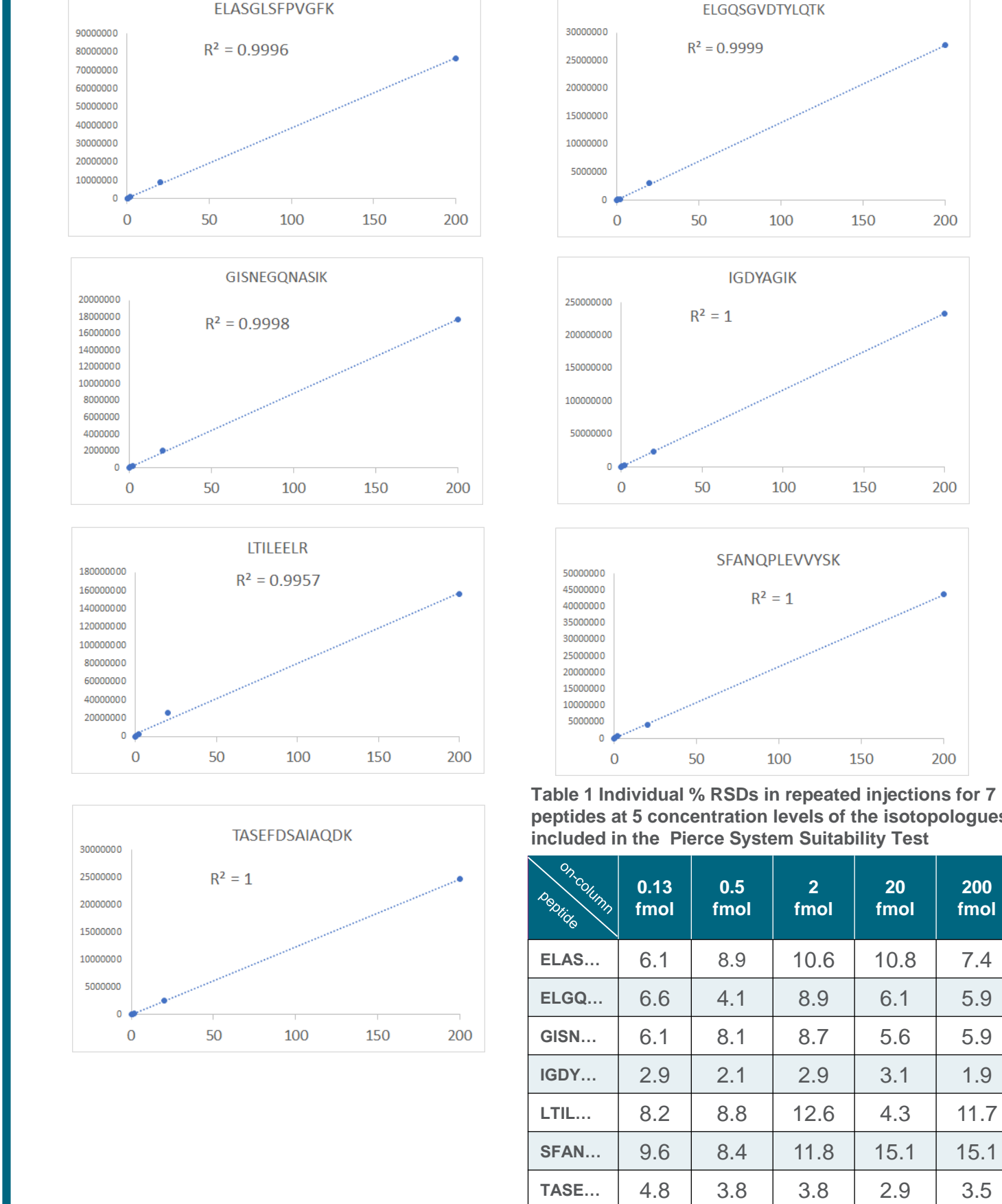


Figure 6 The Pierce LC-MS/MS System Suitability Standard (7 x 5 Mix) enables users to assess performance (sensitivity and dynamic range) of LC-MS/MS systems. The mixture contains 7 peptides, each having 5 isotopologue sequences present in a dilution series. Here the 7x5 Mix was used to evaluate performance of FAIMS Pro with TSQ Altis MS. The mixture was prepared according to Pierce protocol and 0.3ug/ul digested plasma was used as a matrix. The results - obtained with optimized CV values and with 4 SRM transitions for each isotopologue - show linearity for all peptides and low %RSDs for repeated injections.

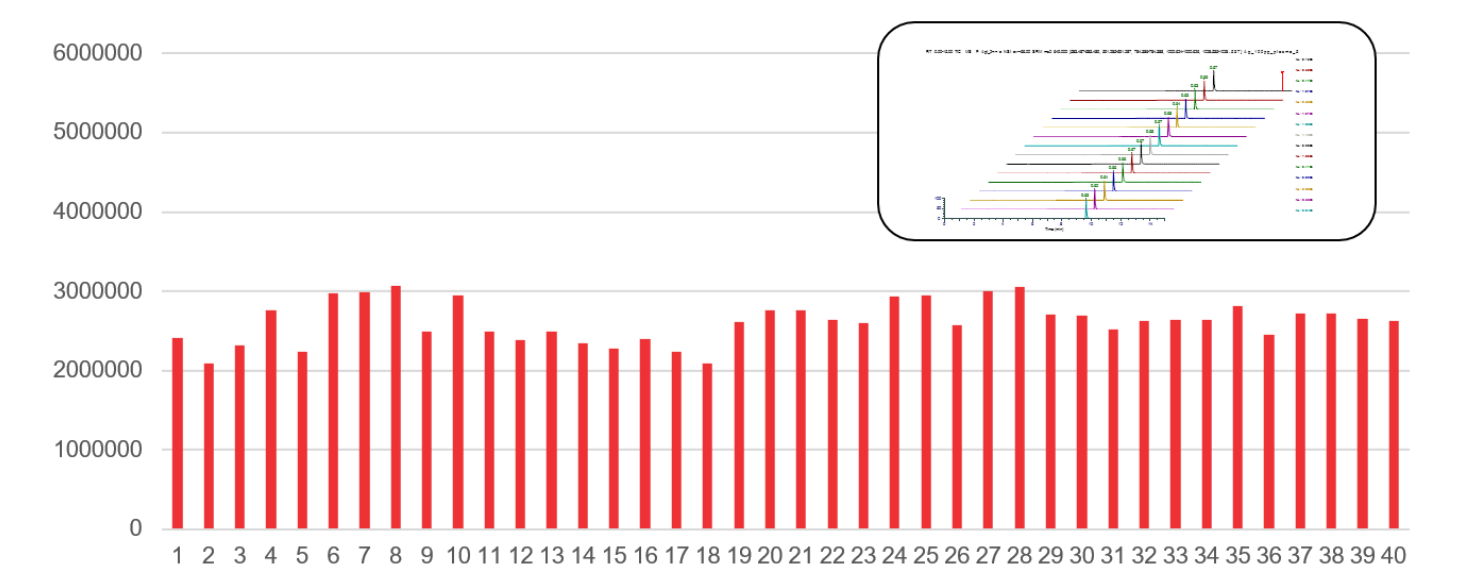


Fig 7 FAIMS Pro – TSQ Altis MS: 40 consecutive injections of Angiotensin in 0.5ug/uL digested plasma. The overall %RSD was <10% across this 2 day experiment (cleaning injections were inserted every 10 analytical injections). The inset show XIC of 15 consecutive injections.

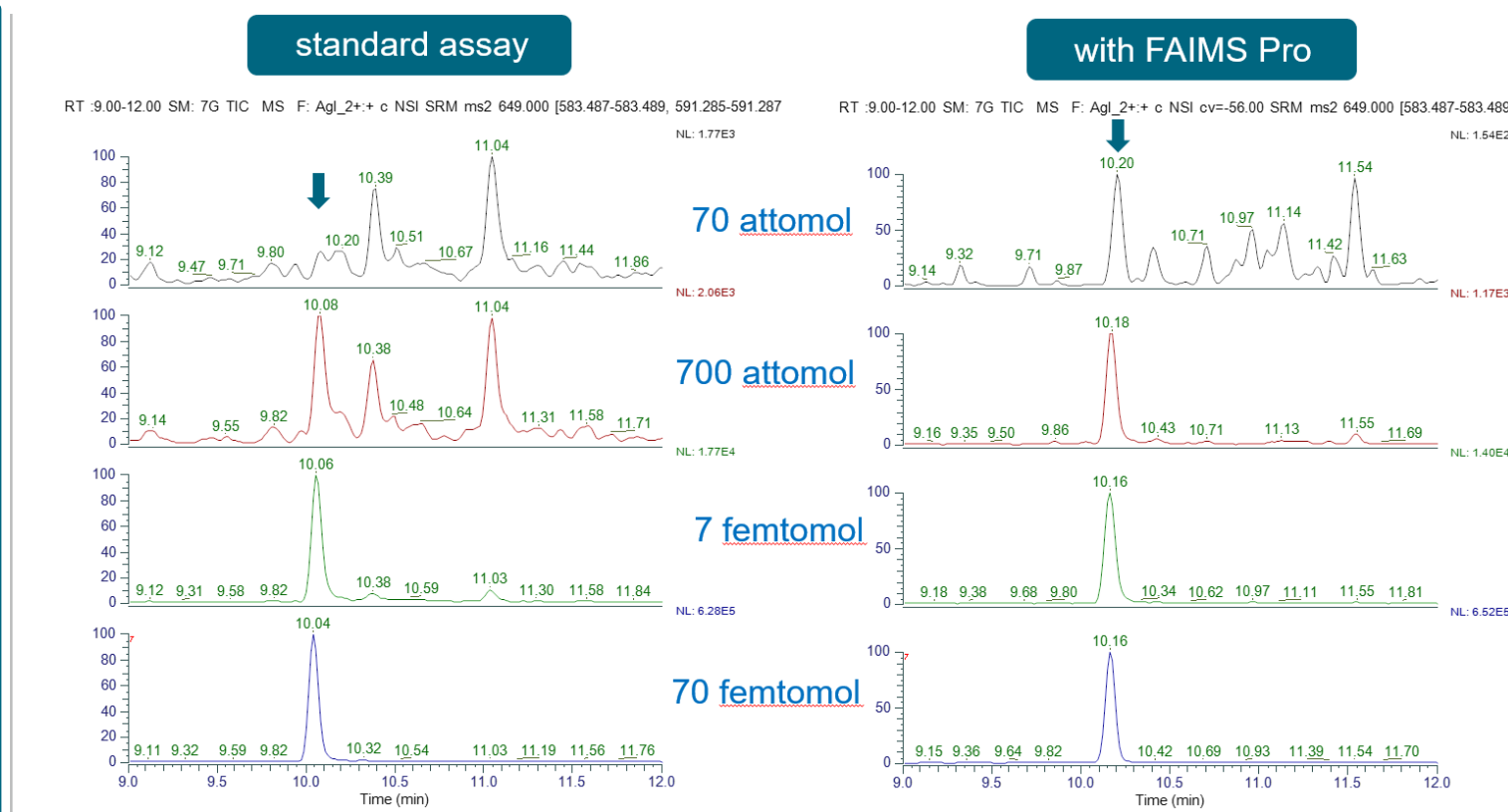


Fig 8 FAIMS Pro – TSQ Altis: Comparison of Angiotensin LC-MS assay with and without FAIMS Pro interface.

CONCLUSIONS

- FAIMS Pro interface coupled to TSQ Altis Triple Quadrupole Mass Spectrometer provides improvement of peptide quantification due to reduced noise and removal of co-eluting interference species.
- FAIMS Pro interface offers orthogonal precursor ion selectivity based on differential gas phase mobility. The Compensation Voltage (CV) setting determines which groups of ions pass the FAIMS Pro interface further to the mass spectrometer. In SRM workflows on triple stage quadrupoles, CV value can be used as another parameter in SRM table to set selective transmission of targeted precursor ion, while suppressing other ions.
- Utilization of FAIMS Pro interface is robust and the analyte signals are reproducible and stable in time as shown by performing consecutive injections over multiple days with low %RSD and by analyzing Pierce System Suitability Standard.

REFERENCES

- Kolakowski BM et al.: Review of applications of high-field asymmetric waveform ion mobility spectrometry (FAIMS) and differential mobility spectrometry (DMS). Analyst. 2007 Sep;132(9):842-64.
- Bonneil E et al. : Enhancement of mass spectrometry performance for proteomic analyses using high-field asymmetric waveform ion mobility spectrometry (FAIMS). J Mass Spectrom. 2015 Nov;50(11):1181-95
- Swearingen KE et al.: High-field asymmetric waveform ion mobility spectrometry for mass spectrometry-based proteomics. Expert Rev Proteomics. 2012 Oct;9(5):505-17.

ACKNOWLEDGEMENTS

Michael MacCoss and Eric Huang of the University of Washington in Seattle and Liz Hecht of Genentech are greatly acknowledged for useful discussions about utilization of the FAIMS Pro technology for the triple quadrupoles peptide quantification workflows. Thermo Fisher Scientific colleagues Scott Peterman, Romain Huguet, Satendra Prasad, Cristina Jacob, Cornelia Boeser, Stephanie Samra and Michael Hudson all kindly provided crucial internal help with this work.

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.



PO65488-EN0419S