

Improving Label-Free Quantification of Plasma and Serum Proteins Using a High-Resolution Hybrid Orbitrap Mass Spectrometer

Maryann S. Vogelsang,¹ Amol Prakash,¹ David Sarracino,¹ Scott Peterman,¹ Bryan Krastins,¹ Jennifer Sutton,¹ Gregory Byram,¹ Gouri Vadali,¹ Shadab Ahmad,¹ Bruno Darbouret,² Mary F. Lopez¹

¹Thermo Fisher Scientific, BRIMS (Biomarker Research Initiatives in Mass Spectrometry), Cambridge, MA, USA;

²Thermo Fisher Scientific, CD Biomarkers, Nimes, France



Overview

Assessing the differences between MS1- and MS2-based label-free relative quantification in a complex plasma matrix using a novel real-time, intelligent acquisition strategy for high-resolution, accurate-mass (HR/AM) global targeted quantification.

Introduction

Label-free mass spectrometry (MS) is an increasingly preferred method for biomarker discovery workflows applied to serum and plasma samples. Given the right conditions, label-free relative quantification is cleaner, simpler, and higher throughput. Resulting differential analysis from these label-free discovery experiments often leads to targeted analyses for verification. High resolution and mass accuracy are critical components to successfully identifying and quantifying peptides in a label-free experiment. Here we present a real-time intelligent acquisition strategy for HR/AM targeted quantification and compare it to relative quantification from MS1 full scan spectra, and introduce a strategy that enables higher confidence in both qualitative and quantitative results in the label-free discovery runs. We propose using HR/AM MS and MS/MS schemes in conjunction with validated spectral libraries for automated method building, data acquisition, verification, and quantification in real-time using novel acquisition schemes.

Methods

Sample Preparation

A protein mixture consisting of eight proteins — cytochrome c (horse), α -lactalbumin (bovine), serum albumin (bovine), carbonic anhydrase (bovine), ovalbumin (chicken), α -S1-casein (bovine), α -S2-casein (bovine), β -casein (bovine) — was prepared at equimolar ratios. The eight non-human proteins were analyzed at 100 fmol on column in a “neat” background as well as 100 fmol on column spiked into a human plasma matrix of 1 μ g on column. The eight proteins were also investigated in the human plasma matrix at varying amounts ranging from 0.5 to 500 fmol each protein on column.

MS Data Acquisition and Analysis

All samples were digested with trypsin and analyzed on a Thermo Scientific™ Q Exactive™ mass spectrometer equipped with a Thermo Scientific™ Nanospray Flex Ion Source. Data was acquired in two steps to simulate traditional workflows. Initial experiments employed unbiased data-dependent MS/MS acquisition resulting in peptide/protein identification as well as building of a spectral library. These initial data-dependent runs were run on both the “neat” conditions of the eight protein mix (without the plasma matrix), and then on a 100 fmol level (each protein) on column in a plasma matrix of (1 μ g plasma on column). These initial data-dependent runs were searched against a modified human database containing the eight additional proteins. The combined results from the discovery experiments were used to build a local spectral library consisting of precursor and product ion m/z values and relative abundance distribution as well as relative retention time values. A highly multiplexed, targeted protein list was created from the spectral library and used for automated data acquisition and processing real-time to facilitate changes to the acquisition scheme. For full description of acquisition method and scheme, please visit poster 131 on Tuesday, by Prakash *et. al.*¹

Thermo Scientific™ Proteome Discoverer™ version 1.3 and Thermo Scientific™ Pinpoint™ version 1.3 software packages were used to analyze both the qualitative and quantitative data. The spectral library resulting from initial runs was used to create a targeted inclusion list and reference information to perform qual/quant determination in real time. Data were acquired and peptide coverage and relative quantification were measured for each of the eight standard proteins. All samples were run in triplicate.

Results

Intelligent Real-Time Data Acquisition

The discovery experiments were performed in an unbiased data-dependent acquisition for the eight protein mixtures in “neat” conditions as well as in a complex plasma matrix. From these initial results, 170 target peptides from the eight proteins were used to build the spectral database, Figure 1. These 170 targets were built into a spectral library look-up table that was used in real-time state modeled acquisition. The look-up table includes the precursor m/z values for the defined charge state as well as the expected retention time window, which are used to initiate product ion spectral acquisition based on the presence of multiple precursor isotopes during the expected elution window (Figure 2).

FIGURE 1. Histogram showing the number of targeted peptides with confident MS2 peak area quantification per femtomolar level of protein mixture. The bars represent the number of confident areas quantified out of the potential 170 targets on the spectral library look-up table.

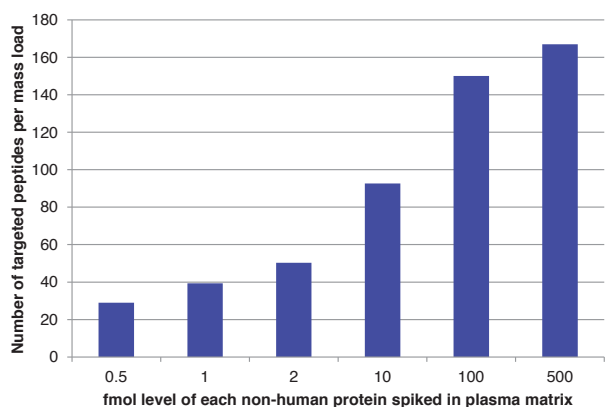
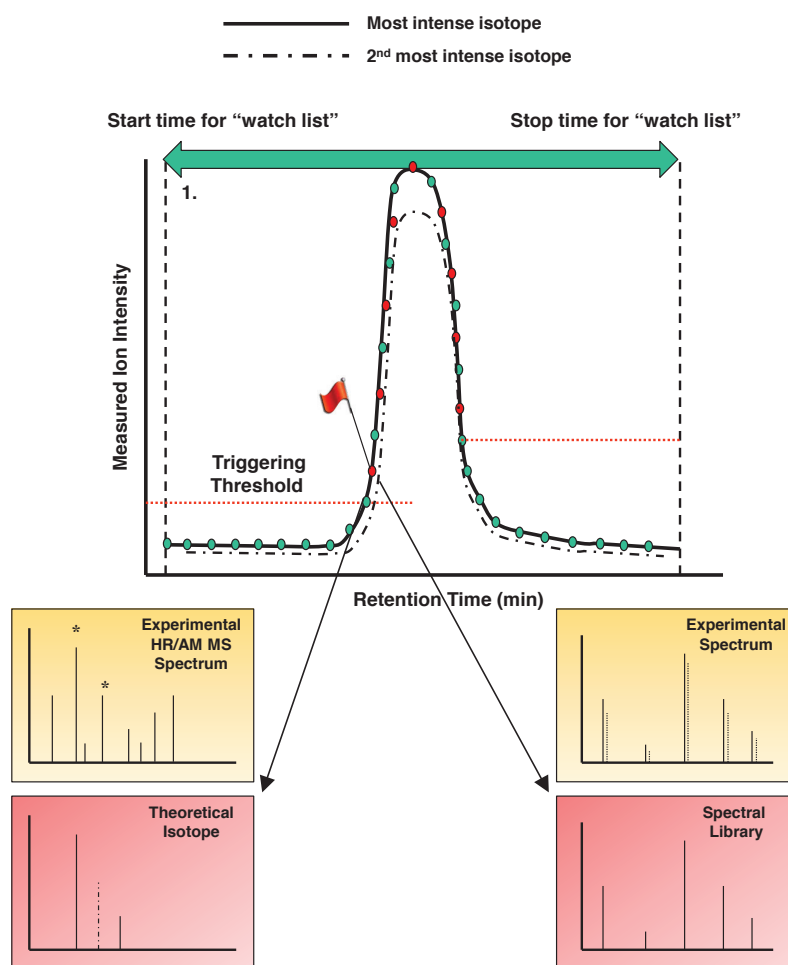


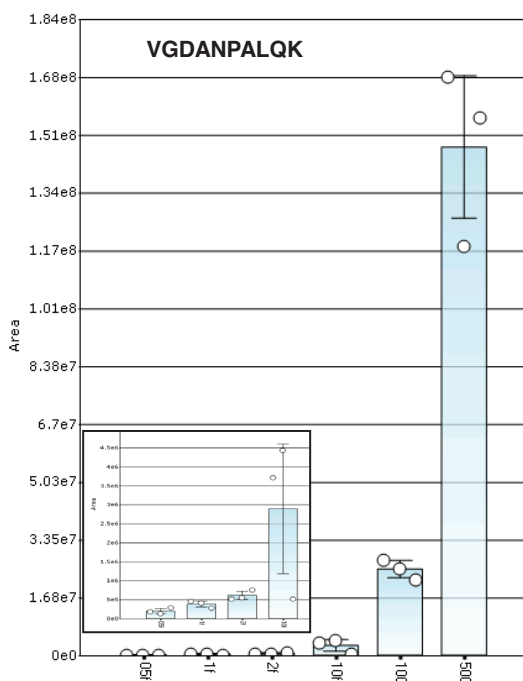
FIGURE 2. Pictorial representation of high IQ data acquisition schemes for targeted peptide quantification using a targeted scanning window, target elution identification, and real-time product ion spectral acquisition. Both precursor and product ion spectral matching is performed to increase the selectivity of data acquisition.



Our current MS-based biomarker discovery studies employ label-free quantification of proteomic data using MS1 extracted ion chromatograms (XIC). In our typical discovery experiments, we get MS1 quantification at signal threshold levels as low as $1e^5$. However, when employing powerful software such as Pinpoint software, we can verify that some of the MS1 quantification may be false positives (demonstrated by loss of multiple isotope peaks or isobaric contaminants).

In this study, we compare the novel real-time state modeled acquisition based MS2 quantification with full scan MS1 (XIC) based quantification (Figure 3). As a general observation, we see that the quality of the MS1 quantification for peptides above 2 femtomoles, the number of targeted peptides we quantify with confidence, is on-par with the MS2-based quantification. However, below 2 femtomoles, the quantity and quality of peptides quantified based on MS1 XIC, drops to nearly 50% of MS2. There is 25–50% false positive quantification at the 0.5–1 fmol level peptides.

FIGURE 3. MS2 Peak area as a function of mass load and their corresponding variance per run for peptide VGDANPALQK. Inset: Expanded view of the 0.5 to 10 fmol level.



Figures 4 and 5 demonstrate the increased confidence that is attained with real-time state-modeled MS2 data quantification. The advantage of this acquisition scheme is evident at levels of proteins and peptides below 10 fmol on column (or at signal thresholds below $1e^4$). There are cases in our study where both MS1 and MS2 do not quantify a species; however, it is important to note that there are significantly fewer false-positives with MS2 quantification.

FIGURE 4. Peak profiles for peptides identified at each femtomolar level, for MS1 quantification (A, C, E) and for MS2 quantification (B, D, F). (A) MS1 peak profiles for peptide VGDANPALQK. (B) MS2 peak profile view of VGDANPALQK. Insets (A) and (B) are expanded views of 0.5 to 2 fmol level. (C) and (D) MS1 and MS2 peak profiles, respectively, for peptide LGEYGFQNALIVR. Insets (C) and (D) are expanded views of 0.5 to 2 fmol level. (E) and (F) MS1 and MS2 peak profiles, respectively, for peptide FALPQYLK. Inset (E) shows an expanded view of probable isobaric contamination from neighboring peak. Inset (F) shows MS2 confidence, 10 fmol level quantification. (G) and (H) MS1 and MS2 peak profiles, respectively, for peptide DM[Oxid]PIQAFLLYQEPVLPVR. Inset (H) shows an expanded view of peak profile for the product ions for the representative peptide at 10 fmol level.

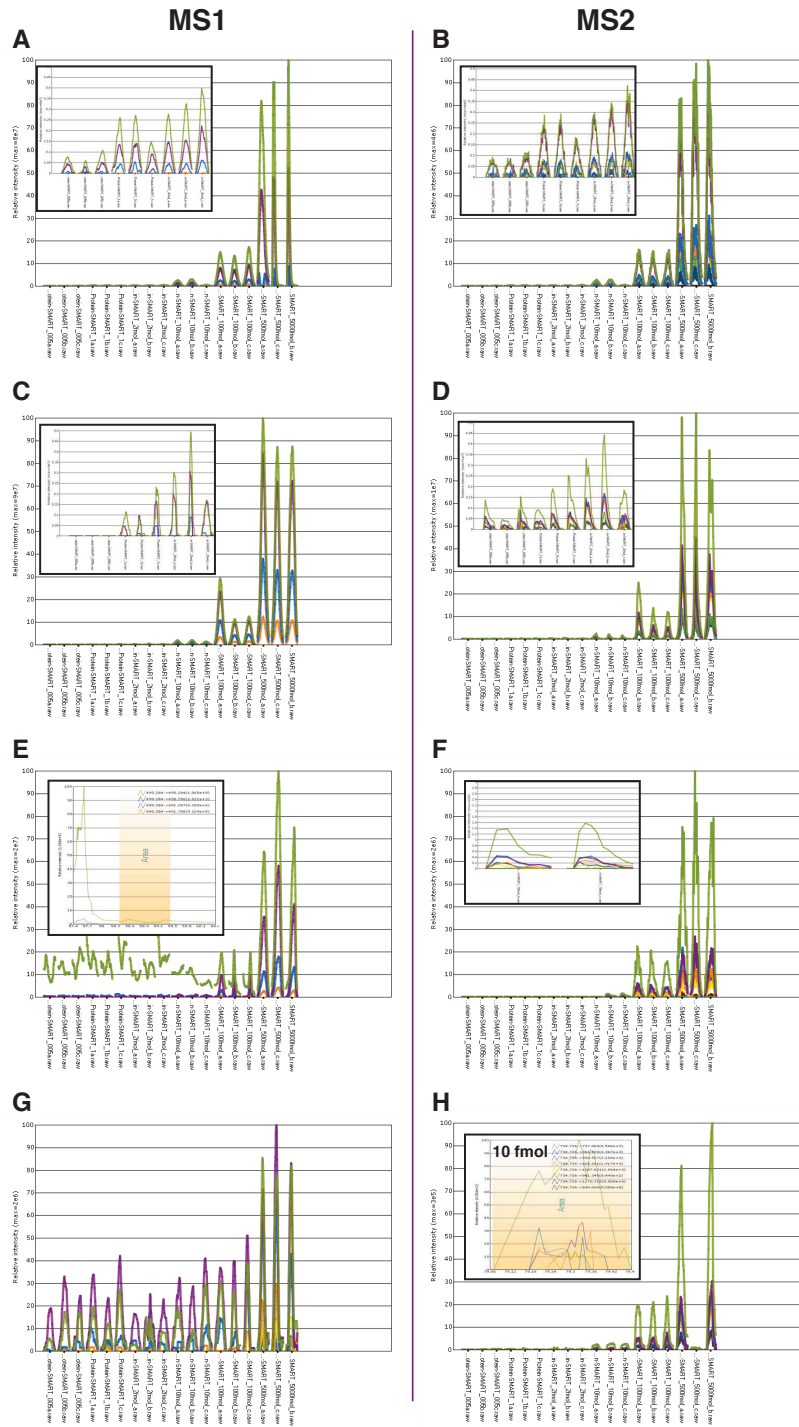
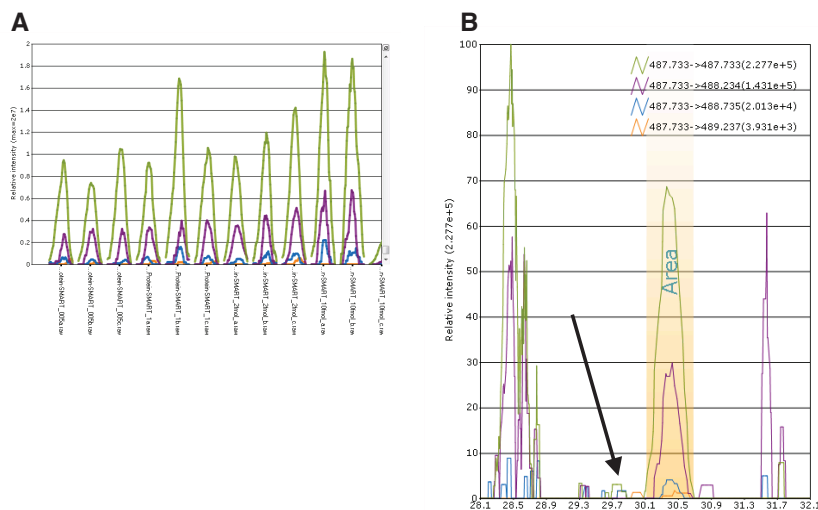


FIGURE 5. (A) False positive MS1 peak profiles for peptide DLGEEHFK at 0.5 to 10 fmol level. (B) XIC for 487.733 isotopes, illustrating probable isobaric contaminants; (arrow) points to observed elution time for peptide DLGEEHFK at higher levels.



Conclusion

- The development of the real-time state modeled data acquisition provides quantification of peptide species at lower concentrations and lower signal thresholds (below $1 e^4$).
- Real-time state modeled data acquisition results in at least 50% more identifications than MS1 quantification below 10 fmol.
- This novel data acquisition scheme provides higher sensitivity and selectivity of peptides in a label-free complex matrix – ideal for a biomarker discovery workflow.

References

1. Prakash, A.; Peterman, S.; Frewen, B.; Kuehn, A.; Ciccimaro, G.; Schroeder, T.; Vasicek, L.; Hood, B.; Bomgardner, R.; Krastins, B.; Sarracino, D.; Byram, G.; Vogelsang, M.; Worboys, J.; Jorgensen, C.; Conrads, T.; Lopez, M. Improving throughput for highly multiplexed targeted quantification methods using novel API-remote instrument control and state-model data acquisition schemes. 61st ASMS Conference on Mass Spectrometry and Allied Topics, Minneapolis, MN, June 9–13, 2013. Poster TP08 – Peptides: Quantitative Analysis I, poster number: 131, Tuesday, Halls B&C.

www.thermoscientific.com

©2013 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific, Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Africa-Other +27 11 570 1840
Australia +61 3 9757 4300
Austria +43 1 333 50 34 0
Belgium +32 53 73 42 41
Canada +1 800 530 8447
China +86 10 8419 3588
Denmark +45 70 23 62 60

Europe-Other +43 1 333 50 34 0
Finland/Norway/Sweden
+46 8 556 468 00
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9434
Italy +39 02 950 591

Japan +81 45 453 9100
Latin America +1 561 688 8700
Middle East +43 1 333 50 34 0
Netherlands +31 76 579 55 55
New Zealand +64 9 980 6700
Russia/CIS +43 1 333 50 34 0
South Africa +27 11 570 1840

Spain +34 914 845 965
Switzerland +41 61 716 77 00
UK +44 1442 233555
USA +1 800 532 4752

Thermo
SCIENTIFIC
Part of Thermo Fisher Scientific