

Maximizing proteome coverage through improved on-line Orbitrap peak determination

Graeme C. McAlister¹, Christian Thoeing², Helene L. Cardasis¹, Alex S. Hebert³, Romain Huguet¹, Philip Remeš¹, Andreas Kuehn², Mike Senko¹, Shannon Eliuk¹, Joshua J. Coon³, Vlad Zabrouskov¹;

¹Thermo Fisher Scientific, San Jose, CA; ²Thermo Fisher Scientific, Bremen, Germany; ³University of Wisconsin, Madison, WI

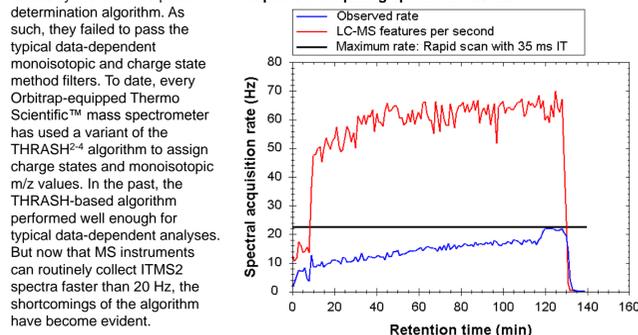
OVERVIEW

Thermo Scientific™ Tribrid™ mass spectrometers are incredibly versatile instruments that combine multiple empowering MS technologies into a single platform. Working together, these technologies can sequence tens of thousands of peptides during a data-dependent LC-MS/MS analysis of a complex peptidic sample. To push sampling depths even further, we have deployed a new peak determination algorithm that identifies hundreds of thousands of additional precursors. To better sample all these additional precursors, we configured the ion trap to collect MS2 spectra at 40 Hz. Together these changes allow us to collect hundreds of thousands of MS2 spectra, which translates into >35% more unique peptide identifications.

INTRODUCTION

The Thermo Scientific™ Orbitrap Fusion Lumos™ Tribrid™ mass spectrometer typically collects ~100,000 ITMS2 spectra during a two hour data-dependent LC-MS/MS analysis of a complex sample (e.g., a tryptic digest of a human lysate). This large pool of MS2 spectra converts to ~30,000 unique peptides. Through this level of proteome coverage is already extensive, exhaustive off-line analysis of the same dataset by HardKlor¹ reveals hundreds of thousands of additional precursors that were never interrogated.

The majority of these unfragmented LC-MS features were never assigned charge states or monoisotopic m/z values. In the past, the THRASH-based algorithm performed well enough for typical data-dependent analyses. But now that MS instruments can routinely collect ITMS2 spectra faster than 20 Hz, the shortcomings of the algorithm have become evident.



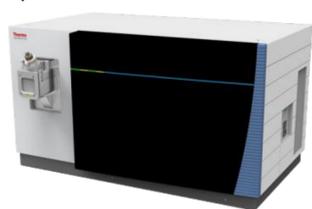
Numerous factors can contribute to the missing or erroneous THRASH-based peak assignments, including overlapping isotopic envelopes and poor ion statistics. Herein we demonstrate the utility of a new on-line Orbitrap peak determination algorithm, which overcomes many of the shortcomings of the old THRASH-based algorithm.

MATERIALS AND METHODS

The Pierce™ HeLa digest protein standard was interrogated using an Orbitrap Fusion Lumos Tribrid mass spectrometer coupled to a Thermo Scientific™ Easy-nLC™ 1000 ultra-high pressure LC. We used a data-dependent method that filtered precursors based upon charge state (2-6), monoisotopic m/z assignment, and dynamic exclusion (20 sec). Unless noted otherwise in the text, ITMS2 spectra were collected at the rapid scan rate, using an automatically determined mass range, and a maximum injection times of 35 ms. The resulting LC-MS/MS data were searched using Thermo Scientific™ Proteome Discoverer™ 2.1 software. The spectra were searched against the UniProt human database, and the peptide spectral matches were filtered to a 1% false-discovery rate using Percolator.

The Thermo Scientific™ Pierce™ Intact Protein Standard mix was also analyzed with an Orbitrap Fusion Lumos Tribrid mass spectrometer, which we coupled to a Thermo Scientific™ Dionex™ UltiMate™ 3000 ultra-high pressure LC operating at 200µl/min. We analyzed the sample with a data-dependent method that consisted of a 15k resolution (@200m/z) Orbitrap MS1 scan followed by data-dependent Orbitrap MS2 scans with the precursor charge state filter set to ≥7.

Figure 2. All the data were collected using an Orbitrap Fusion Lumos Tribrid mass spectrometer.



RESULTS

Initial characterization of Advanced Peak Determination algorithm

The Advanced Peak Determination (APD) algorithm boasts a suite of new features. These include the ability to annotate overlapping isotopic envelopes, improvements to the pattern matching filters used to assign the charge states and monoisotopic m/z values (e.g., the Patterson filter and the average model correlation), and a function that correlates assignments across the entire charge envelope of a given precursor (i.e., charge state deconvolution).

Between back-to-back LC-MS/MS analyses we alternated between the standard THRASH-based algorithm and the new APD algorithm.

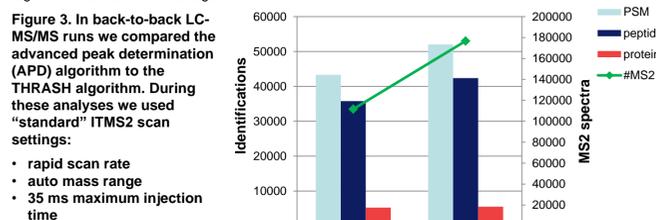


Figure 3. In back-to-back LC-MS/MS runs we compared the advanced peak determination (APD) algorithm to the THRASH algorithm. During these analyses we used “standard” ITMS2 scan settings:

- rapid scan rate
- auto mass range
- 35 ms maximum injection time

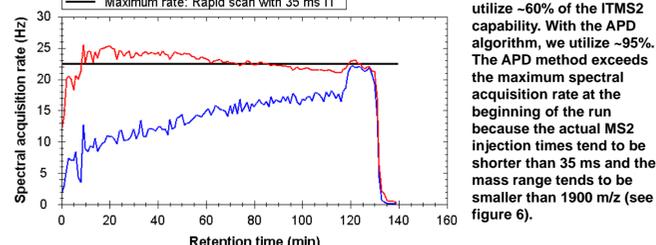


Figure 4. With the THRASH-based algorithm, we only utilize ~60% of the ITMS2 capability. With the APD algorithm, we utilize ~95%. The APD method exceeds the maximum spectral acquisition rate at the beginning of the run because the actual MS2 injection times tend to be shorter than 35 ms and the mass range tends to be smaller than 1900 m/z (see figure 6).

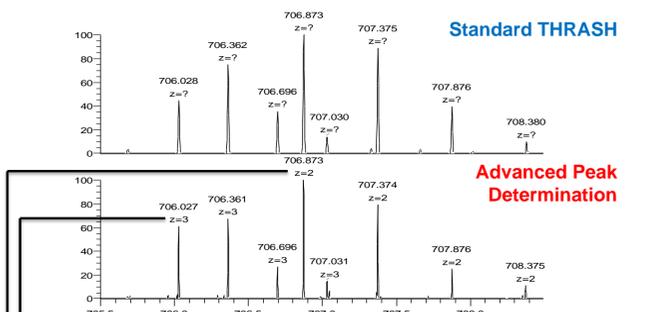


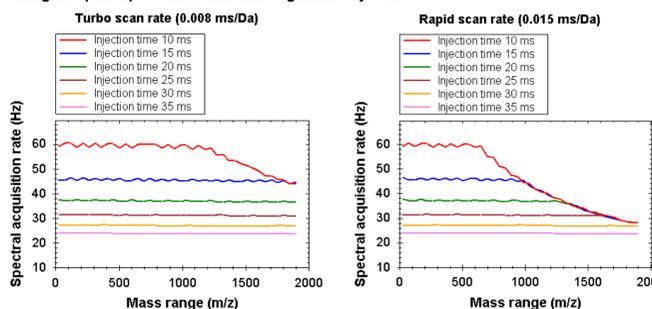
Figure 5. An example FTMS1 mass range from the LC-MS/MS experiments where APD was disabled (top), and where APD was enabled (bottom). The mass range contains overlapping isotopic envelopes that only the APD algorithm can accurately identify.

	Monoisotopic m/z	Charge	Sequence	XCorr
→ Precursor #1	706.027	3	ASLLQNSTNEQLQIHYK	3.568
→ Precursor #2	706.873	2	EPALNEANLSNLK	3.385

Optimization of the ITMS2 scan settings

When the pool of available precursors was limited, it made sense to use the excess MS cycle time to collect higher quality MS/MS spectra at a slower acquisition rate. Now that we have a much larger population of precursors to interrogate, we can afford to collect more MS2 spectra at a faster rate.

Figure 6. The ITMS2 spectrum acquisition rate is determined by the analyzer scan rate, maximum injection time, and scan range. On the Tribrid MS, ion injection occurs concurrently with m/z analysis of the preceding scan. As such, these parameters work together to determine the maximum spectrum acquisition rate. For these graphs, ions were isolated using the quadrupole mass filter and fragmented by HCD.



An optimal LC-MS/MS method needs to weigh the counterbalanced goals of spectral quality and spectral acquisition rate. As ion injection time decreases, the MS2 spectral acquisition rate increases but the ion statistics in the resulting MS2 spectra decrease. Over the course of many LC-MS/MS analyses, we optimized various instrument parameters, including ITMS2 maximum injection time, MS2 mass range, MS2 scan rate, FTMS1 resolution, and MS cycle time.

Figure 7. Varying the maximum ITMS2 injection during a 2 hour LC-MS/MS analysis. The scan rate was turbo and the mass range was 200-1400 m/z.

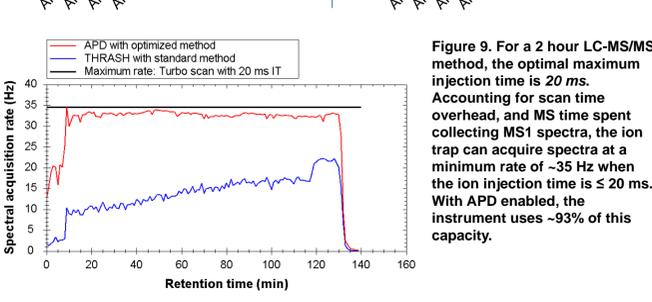
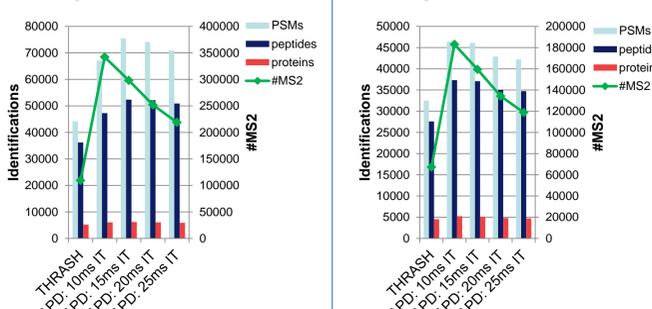


Figure 8. Varying the maximum ITMS2 injection during a 1 hour LC-MS/MS analysis. The scan rate was turbo and the mass range was 200-1400 m/z.

Figure 9. For a 2 hour LC-MS/MS method, the optimal maximum injection time is 20 ms. Accounting for scan time overhead, and MS time spent collecting MS1 spectra, the ion trap can acquire spectra at a minimum rate of ~35 Hz when the ion injection time is ≤ 20 ms. With APD enabled, the instrument uses ~93% of this capacity.

Replicate analyses using APD LC-MS/MS

For LC-MS/MS methods with APD enabled, we found that the following ITMS2 settings nicely balanced sensitivity and robustness: 20 ms maximum injection time, rapid scan rate, and an automatic mass range. Using these settings, we performed replicate (n=3) LC-MS/MS analyses comparing APD on vs. APD off. For both conditions, we injected 1 µg of a tryptic HeLa digest, and we performed both 1 and 2 hour LC gradients. During the THRASH-based LC-MS/MS method we used our “standard” ITMS2 scan settings: 35 ms maximum injection time, rapid scan rate, and an automatic mass range.

Figure 10. During a 2 hr LC-MS/MS method we identified >35% more unique peptides with APD. Or as an alternative, with APD enabled we identified an equivalent number of unique peptides with an LC-MS/MS method that is half as long as the 2 hr THRASH-based LC-MS/MS method.

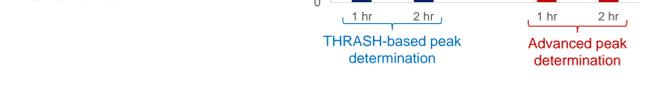


Figure 10. During a 2 hr LC-MS/MS method we identified >35% more unique peptides with APD. Or as an alternative, with APD enabled we identified an equivalent number of unique peptides with an LC-MS/MS method that is half as long as the 2 hr THRASH-based LC-MS/MS method.

Figure 11. We compared the overlap between 3 x APD 2 hr LC-MS/MS analyses to the overlap between 3 x THRASH-based 2 hr LC-MS/MS analyses. The number of unique peptides observed in all three replicates increased by ~50% with APD (26,158 vs. 38,045).

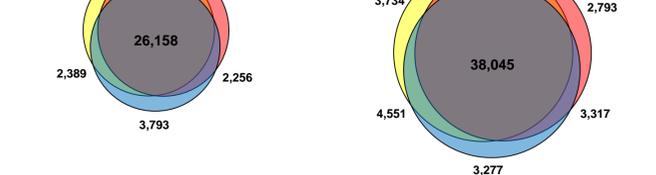


Figure 12. The different populations in the Venn diagrams above (peptides seen in 1, 2, or all 3 replicates) were distributed on a histogram by precursor intensity. In this figure the three populations are stacked (i.e., replicate 1 = replicate 1, replicate 2 = 1+2, and replicate 3 = 1+2+3), such that the entire histogram represents all the peptides identified in all the replicates.

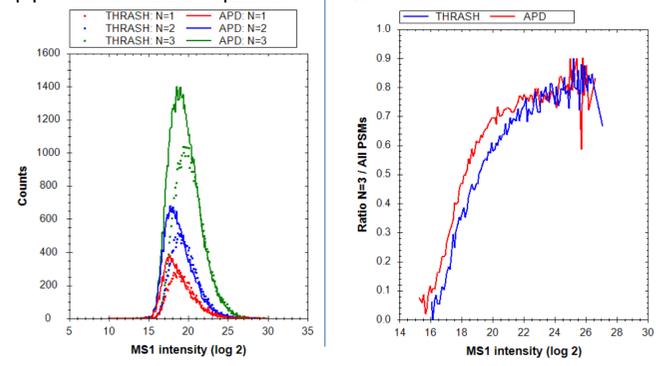


Figure 12. The different populations in the Venn diagrams above (peptides seen in 1, 2, or all 3 replicates) were distributed on a histogram by precursor intensity. In this figure the three populations are stacked (i.e., replicate 1 = replicate 1, replicate 2 = 1+2, and replicate 3 = 1+2+3), such that the entire histogram represents all the peptides identified in all the replicates.

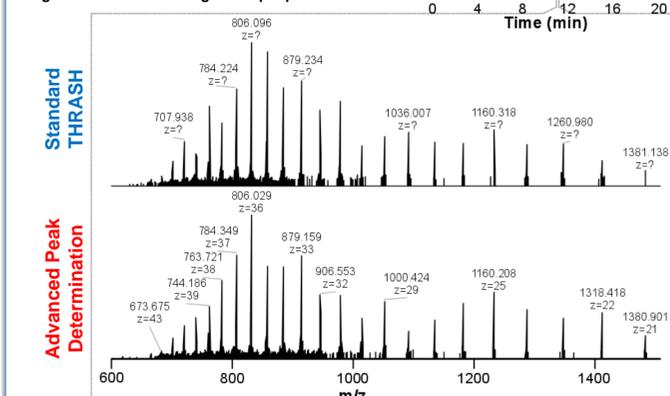
Figure 13. Using the distributions in figure 12, at each intensity bin we calculated the ratio between the number of peptides identified in all three replicates over the total number of peptides. Based upon this analysis, the limit of reproducible detection (where > 50% of all the peptides identified were observed in all three replicates) is ~2-3x lower with APD.

TOP-DOWN RESULTS

Demonstration of the utility of APD for top-down analysis

The APD algorithm's ability to identify overlapping isotopic envelopes is the main force behind most of the gains we observed during the LC-MS/MS analyses of peptide samples. However, the APD algorithm also has improved charge state assignment functions, including the ability to correlate assignments across the entire precursor charge envelope. These other APD advancements greatly improve large biomolecule charge state assignment.

Figure 14. We analyzed the Pierce intact protein standard mix by LC-MS/MS. FTMS1 spectra were collected at 15k and were the summation of 5 scans. In back-to-back runs we tested APD on and off. With APD on, the instrument can accurately identify highly charged and complex protein envelopes. Thanks to the improved charge state assignment abilities of the APD algorithm, we can execute more advanced top-down methods – such as, MS/MS fragmentation of one charge state per precursor.



CONCLUSIONS

- The advanced peak determination (APD) algorithm identifies hundreds of thousands of additional precursors in Orbitrap spectra for data-dependent analysis.
- We observe a large increase in the unique peptide identifications when we configure the quadrupole ion trap to favor a faster MS/MS rate that more effectively samples all the additional precursors.
- With APD and the optimized ITMS2 method, we collect ~250,000 MS2 spectra during a 2 hr LC-MS/MS method. This converts into >45,000 unique peptide IDs, which is a >35% improvement over the conventional THRASH-based approach.
- This improved sample coverage translates into better run-to-run reproducibility. Such that, the limit of reproducible identification is ~2-3x lower with APD.
- APD also enables improved large molecule charge state assignment and in turn top-down data-dependent decisions.

REFERENCES

1. Hoopmann MR, Finney GL, MacCoss MJ. High speed data reduction, feature selection, and MS/MS spectrum quality assessment of shotgun proteomics datasets using high resolution mass spectrometry. *Anal. Chem.* 2007, 79, 5630-5632.
2. Senko MW, Beu SC, McLafferty FW. Automated assignment of charge states from resolved isotopic peaks for multiply charged ions. *J. Am. Soc. Mass Spectrom.* 1995, 6, 52-56.
3. Senko MW, Beu SC, McLafferty FW. Determination of monoisotopic masses and ion populations for large biomolecules from resolved isotopic distributions. *J. Am. Soc. Mass Spectrom.* 1995, 6, 229-233.
4. Horn DM, Zubarev RA, McLafferty FW. Automated reduction and interpretation of high resolution electrospray mass spectra of large molecules. *J. Am. Soc. Mass Spectrom.* 2000, 11, 320-332.

ACKNOWLEDGEMENTS

The authors would like to thank Michael Westphal, Jae Schwartz, John Syka, Jesse Canterbury, Chris Mullen, and Rosa Viner for helpful conversations.

© 2017 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.