

# Performance metrics and example applications: Orbitrap mass spectrometry at resolving power 1M

Jesse Canterbury<sup>1</sup>, Michael Senko<sup>1</sup>, Seema Sharma<sup>1</sup>, Tim Stratton<sup>1</sup>, Vlad Zabrouskov<sup>1</sup>, Eduard Denisov<sup>2</sup>, Alexander Makarov<sup>2</sup> Thermo Fisher Scientific, <sup>1</sup>San Jose, CA, USA and <sup>2</sup>Bremen, Germany

## ABSTRACT

**Purpose:** This work surveys key performance problems and their solutions in Orbitrap mass spectrometry at high resolving powers and presents example applications benefiting from high-resolution Orbitrap analysis.

**Methods:** A Thermo Scientific™ Orbitrap Fusion Lumos™ Lumos™ Tribrid™ mass spectrometer was modified by (1) ultra-precisely machined Orbitrap electrodes; and (2) carefully controlled injection conditions.

**Results:** Improvements in performance enabling routine analysis at resolving power 1,000,000 (at  $m/z$  200), are presented, along with results from applications in metabolomics and lipid fluxomics.

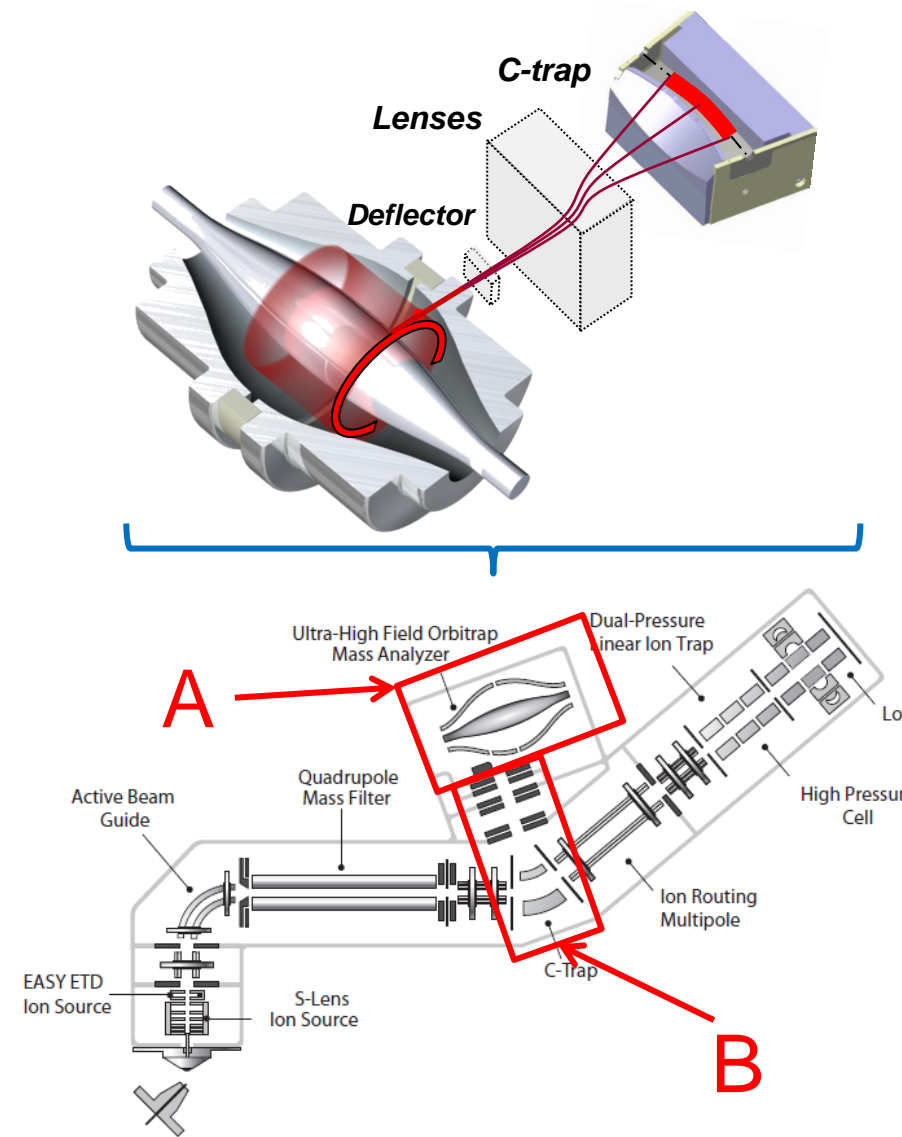
## INTRODUCTION

Since its commercial introduction in 2005, the Orbitrap mass analyzer has been widely adopted by the mass spectrometry community. Historically, applications in which Orbitrap analysis has been particularly useful are those which require high resolving power. For example, in many proteomics studies, resolution of 60k or 120k (at  $m/z$  200) is more than sufficient to accurately determine the monoisotopic mass of multiply charged peptides across the mass range of interest. However, some applications require higher resolutions and until now they remained served only by Fourier Transform Ion Cyclotron Resonance (FT-ICR) instruments. Here we describe an Orbitrap-based mass spectrometer capable of dependable and routine analysis at a resolving power of 1M. Performance criteria are discussed and applications examples are provided.

## MATERIALS AND METHODS

**Samples.** For calibration and characterization, a standard calibration solution was used, which consisted of n-butylamine, caffeine, MRFA, and Ultramark 1621 dissolved in 50:49:1 acetonitrile, methanol, and formic acid. For the metabolite standards experiment, 24 common metabolites were mixed at a concentration of 100 pg/μL. 20 μL were then injected on to a Thermo Scientific™ Hypersil GOLD™ column (100x2.1 mm). Analytes were separated using a 20-min gradient starting at 5% buffer B (99.9:0.1 acetonitrile:formic acid) and increasing to 35% B (buffer A was 99.9:0.1 water:formic acid), at a flow rate of 300 μL/min. For the flavonoid and lipid fluxomics work, sample information can be found in the referenced presentations.

**Mass spectrometry.** All mass spectrometry was performed on a Orbitrap Fusion Lumos Tribrid mass spectrometer equipped with the 1M Option (Fig. 1). The 1M Option consists of ultra-precisely machined high-field Orbitrap electrodes used in combination with a C-trap and injection optics with improved manufacturing tolerances. The UHV system was baked at a temperature of approximately 120°C until the base pressure was  $\leq 5 \times 10^{-11}$  torr. Applied voltages were calibrated according to an automated calibration routine as described below.

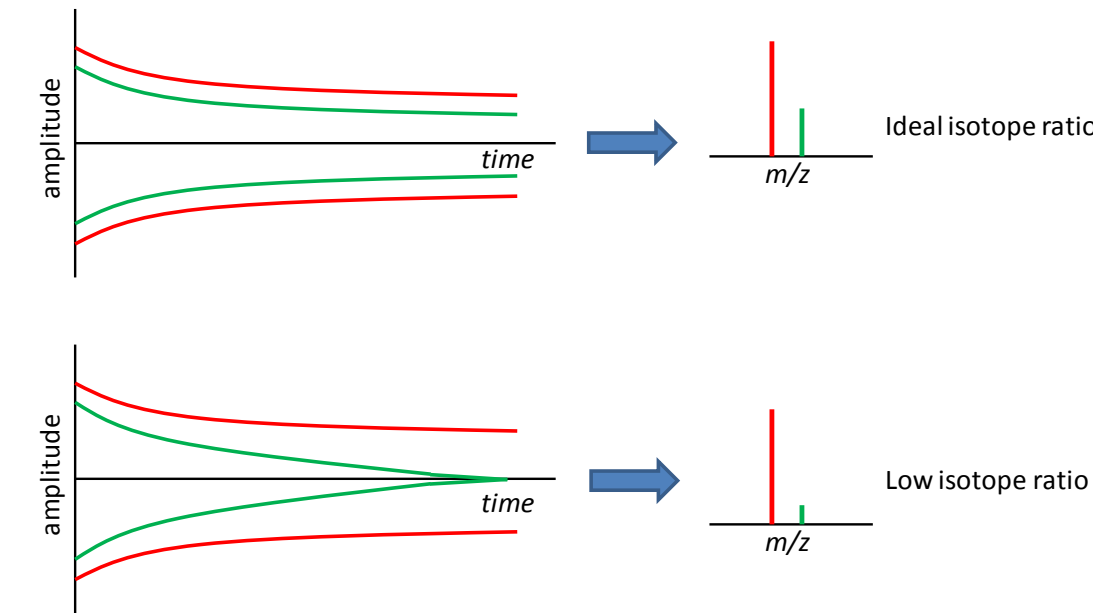


**FIGURE 1.** Orbitrap Fusion Lumos mass spectrometer equipped with the 1M Option: (A) high-precision Orbitrap electrodes; (B) C-trap and injection optics with improved manufacturing tolerances.

## RESULTS

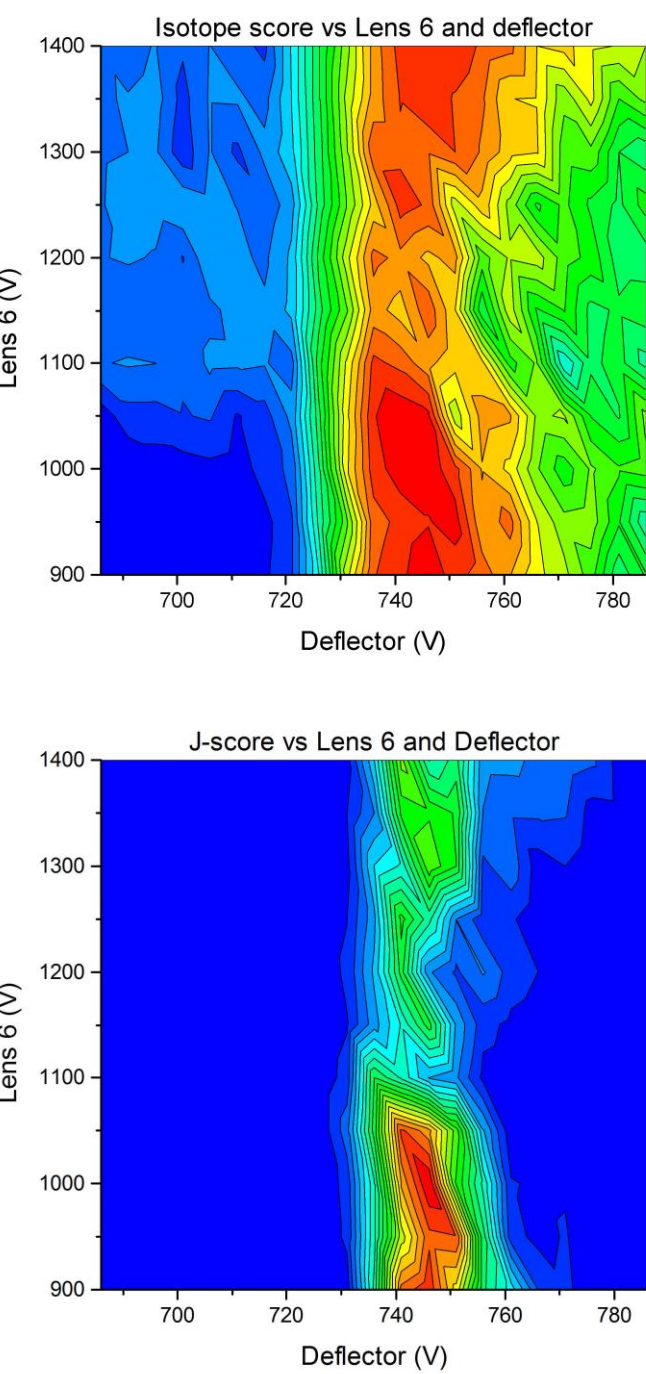
### Performance considerations and validation

In FTMS measurements, ion cloud coherence determines the length of time over which useful measurements can be made. Cloud coherence can be compromised especially for less dense ion clouds, and this can lead to intensity-dependent decay. This in turn can lead to inaccurate isotope abundances, as shown below.



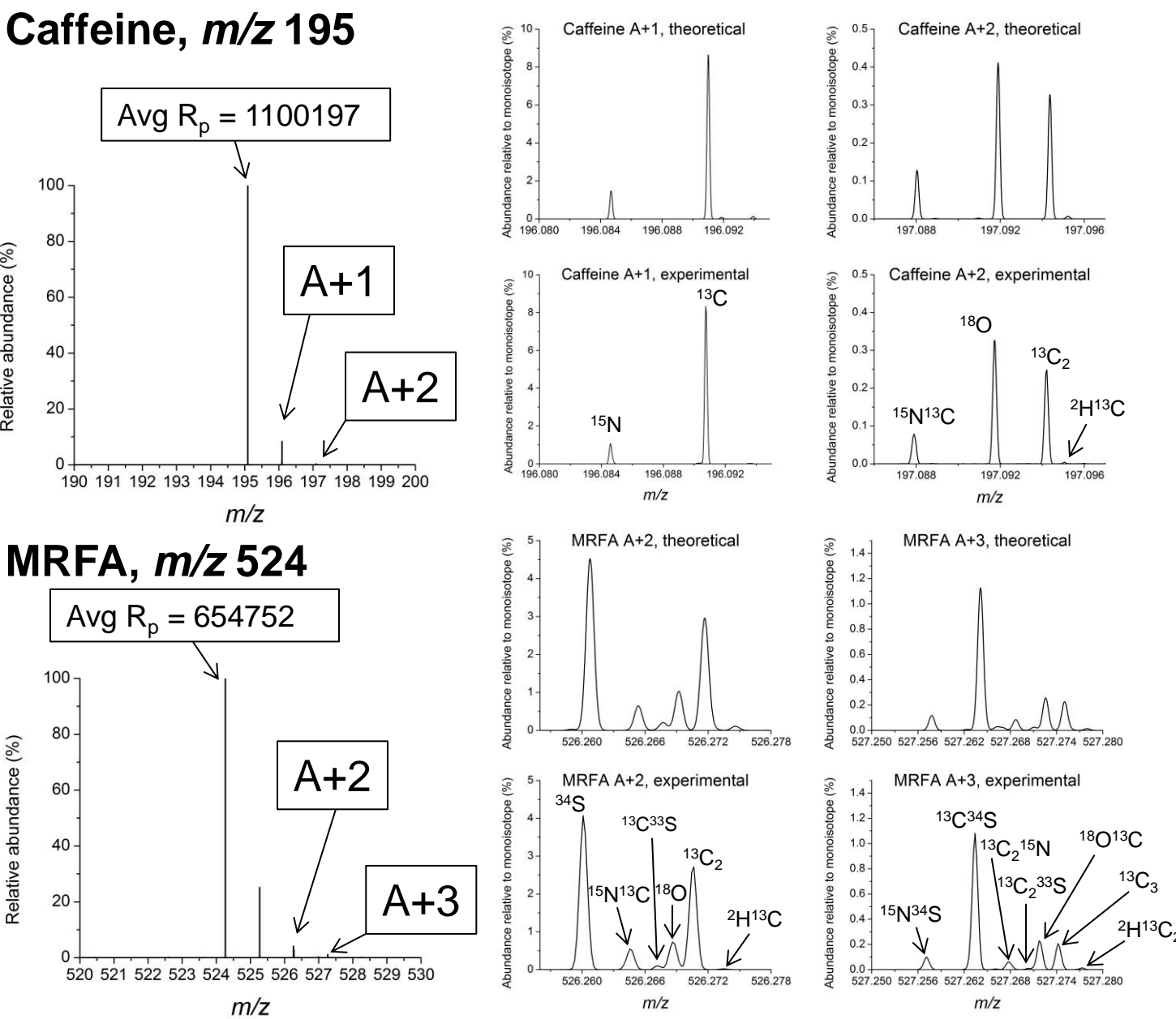
**FIGURE 2.** Intensity-dependent signal decay can lead to inaccurate isotopic abundances for smaller ion clouds (green) relative to larger (red).

The problem described above is worse for longer transients, i.e., high resolving powers. To ensure quantitative accuracy and adherence to all performance objectives, we employ an automated calibration routine. This builds upon the routine previously described [1], which employs a genetic algorithm and tailored scoring in order to navigate the multidimensional optimization surface that arises in C-trap/Orbitrap calibration.

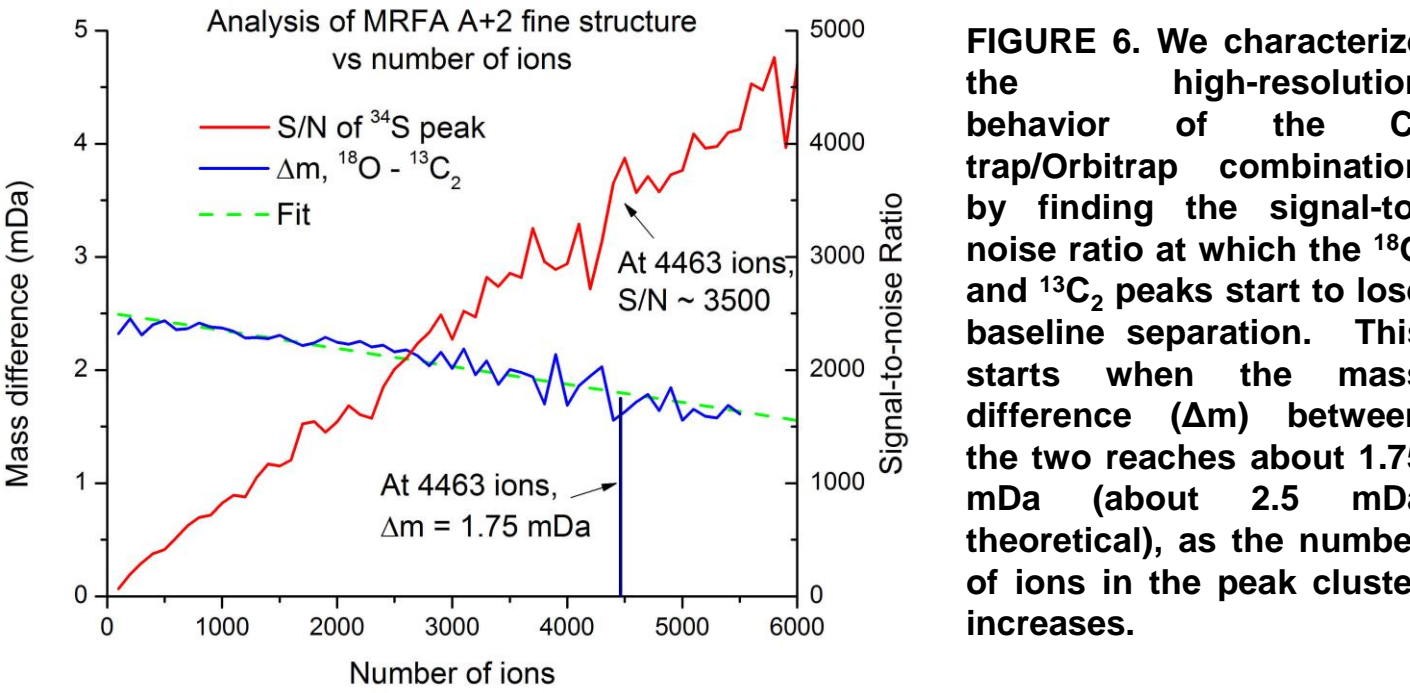
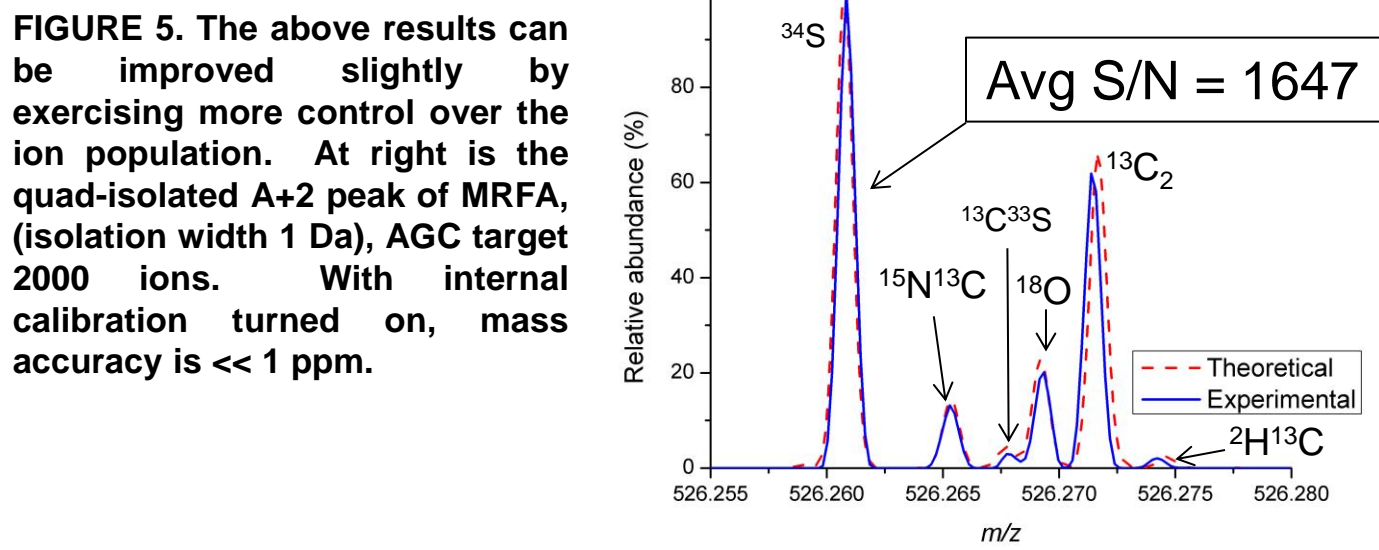


**FIGURE 3.** Top left, complex surface representing the response of an isotope score (a composite of isotope fidelity measurements for caffeine, MRFA, and UM1522) to changes in two important optics elements, the deflector and an adjacent lens. While complex, this 2-D surface captures only a portion of the complexity of the entire multidimensional surface, which includes responses from 6 other optics elements.

Bottom left, this complexity can be significantly reduced by using a more selective score (see ref [1]), which makes the true optimum more clear.



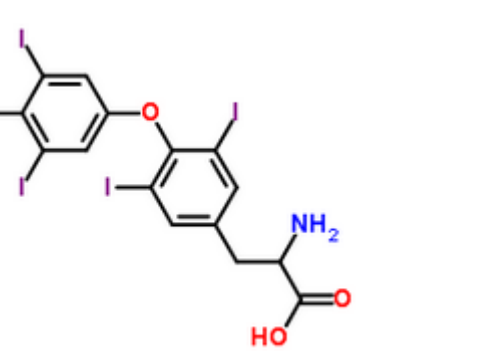
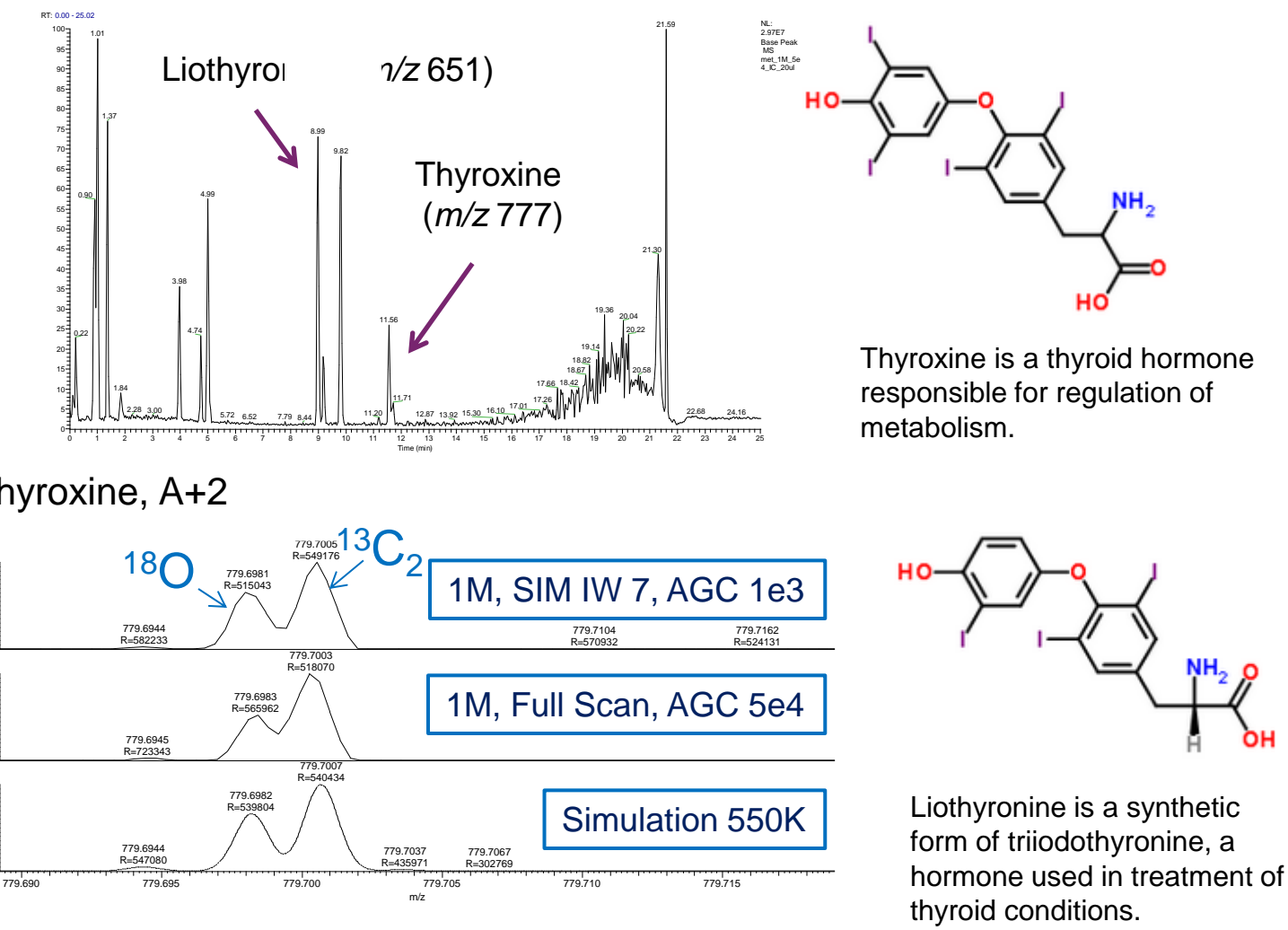
**FIGURE 4.** With the improvements shown described in methods and with proper calibration, good results can be obtained as shown above. Each species is quad-isolated (20  $m/z$  isolation width), AGC target 5e4 ions.



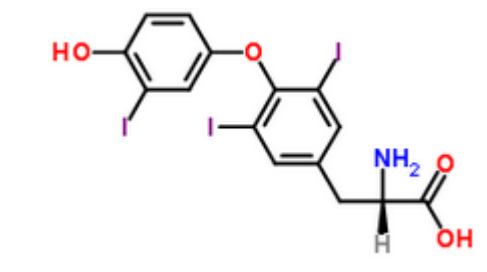
## Example applications

The figures below show fine isotopic structure revealed in a metabolites standard, on an LC timescale.

### LC/MS of metabolite standards



Thyroxine is a thyroid hormone responsible for regulation of metabolism.

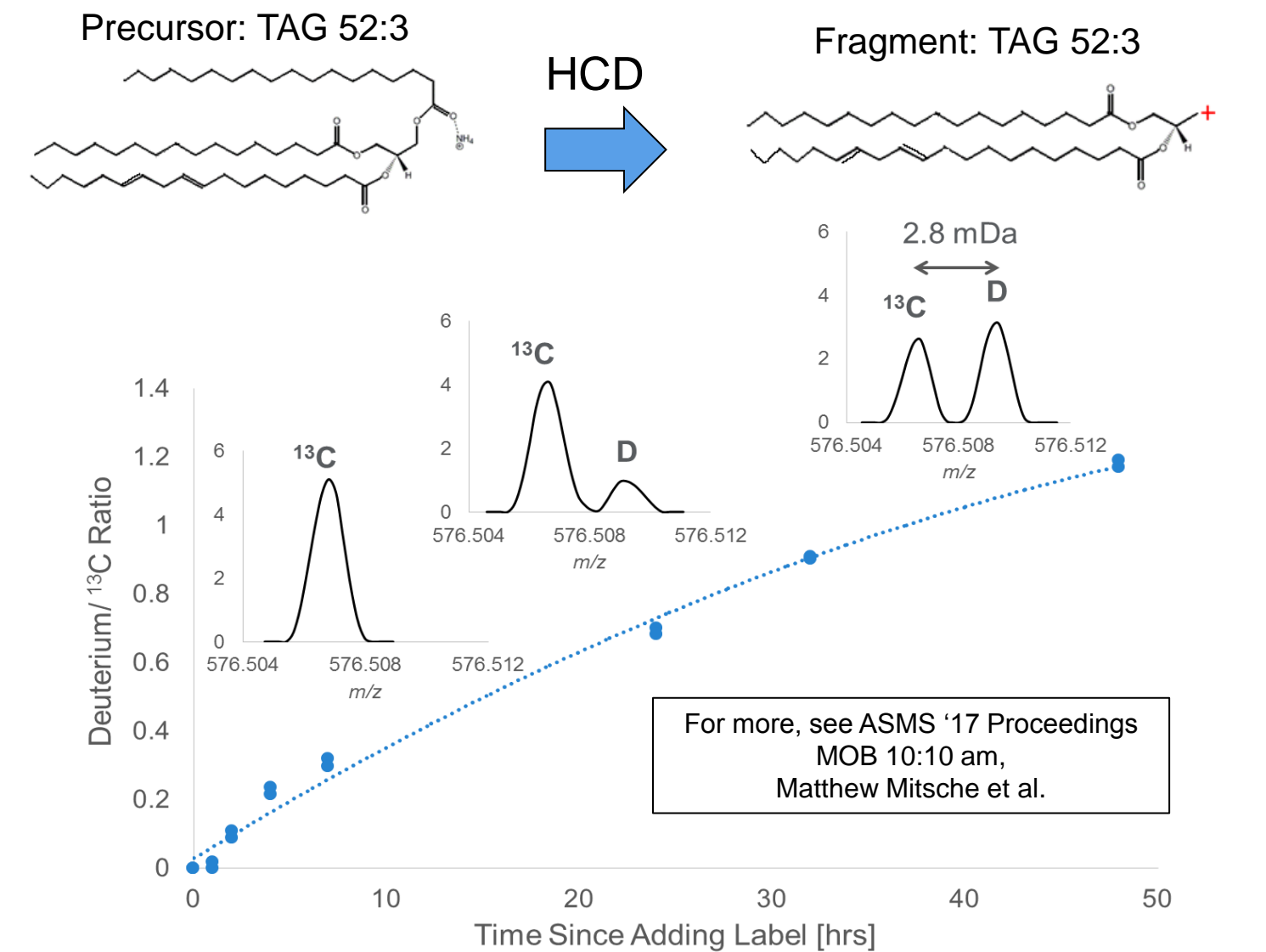


Liothyronine is a synthetic form of triiodothyronine, a hormone used in treatment of thyroid conditions.

### Lipid Fluxomics

In *fluxomics*, the objective is the quantitation of functional molecular turnover over many different molecular species of interest.

The figure below describes measurements of lipid biosynthesis. Normally such measurements are made using tritium labeling; however, very few laboratories are able to practice this method due to strict controls on radioactive materials and handling. Here deuterium labeling is used, and the labeled and unlabeled forms can be baseline separated using ultra-high resolution Orbitrap MS.



**FIGURE 9.** Intact deuterium-label triacylglyceride (TAG 52:3) is fragmented to yield a diacylglyceride (DAG 34:1), which retains a double-bond. The fragment DAG is at lower  $m/z$  where higher resolving power allows the <sup>13</sup>C and D peaks to be easily separated.

## CONCLUSIONS

- Using ultra-precisely machined Orbitrap electrodes and employing exact control over ion injection conditions, transients of 2 seconds yield good performance at a mass resolving power of 1,000,000 (at  $m/z$  200).
- Optimized calibration and tuning provide accurate isotope fidelity and robust space charge performance
- Data obtained with the 1M Option provide important information about a variety of research topics.

## REFERENCES

- J. Canterbury et al., Poster M309, ASMS 2013, "Automatic Isotope Ratio Optimization on High-field Orbitrap Mass Analyzers."

## ACKNOWLEDGEMENTS

We are grateful to Matthew Mitsche of the University of Texas Southwestern Medical Center for examples from his lipid fluxomics work.

We gratefully acknowledge Dae-Eun Lee for valuable assistance during the course of this work. We thank Hartmut Kuipers and Oliver Lange for help with data processing. Finally we thank Raman Mathur and Chad Weisbrod for helpful discussions.

## TRADEMARKS/LICENSEING

© 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.