Developments in Orbitrap mass spectrometry on a modified Tribrid mass spectrometer

Jesse Canterbury¹, Graeme McAlister¹, Michael Senko¹, Romain Huguet¹, Aaron Robitaille¹, Arne Kreutzmann², Konstantin Aizikov², Dmitry Grinfeld², and Alexander Makarov² ¹Thermo Fisher Scientific, San Jose, California, USA; ²Thermo Fisher Scientific, Bremen, Germany

ABSTRACT

Purpose: Improved instrument performance at elevated pressures and for higher m/z analytes; improved speed of analysis for TMT workflows

Methods: We modified a standard Thermo Scientific[™] Orbitrap[™] Fusion Lumos[™] Tribrid[™] mass spectrometer. Modifications included software and hardware changes, as described below.

Results: (1) Sensitivity for larger analytes is improved by raising the pressure in the ion routing multipole (IRM) and slightly modifying transmission characteristics. (2) Resolution or speed for TMT work can be improved by employing TurboTMT, which uses the ΦSDM algorithm, delivering higher resolution than normal FT methods for the same transient length.

INTRODUCTION

Improving the acquired information per unit time is an important goal for any LC-MS workflow. Information flux is improved as the breadth and quality of data collected on a single instrument expands. In this work, we describe modifications to a Tribrid system for the purpose of extending the overall utility of the system in a variety of applications, from shotgun proteomics to large molecule analysis. These modifications include increasing the mass range, enabling a wider range of system pressures, and deploying new spectral annotation algorithms.

MATERIALS AND METHODS

Sample Preparation. In most cases, characterization was done using Thermo Scientific[™] Pierce[™] FlexMix[™] calibration solution. For the protein measurements in Figure 4, carbonic anhydrase from bovine erythrocytes (Sigma-Aldrich, St. Louis) was dissolved in 50:49.9:0.1 water:methanol:formic acid to a concentration of 1 µM. TMT work was done with the Thermo Scientific[™] Pierce[™] TMT11plex yeast digest standard.

Instrument Modifications. We modified a standard Orbitrap Fusion Lumos Tribrid mass spectrometer as described below, and shown in Figure 1. Hardware modifications included elevating the pressure in the IRM as well as Enhanced Vacuum Technology, a proprietary modification of the vacuum system. Software modifications included integration of the Φ SDM algorithm [1] into the instrument control software.



Figure 1. Modified Orbitrap Fusion Lumos Tribrid mass spectrometer. Changes discussed in this work are in the IRM and C-trap/Orbitrap area.

RESULTS

Operation at higher IRM pressure

the cell.

0.020

Figure 2. As shown at right, IRM pressure can be accurately regulated across a wide range of pressures, from 1 mtorr up to 20 mtorr.

0.015 -

5.6x10^{-1*}

5.4x10^{-1*}

Figure 3. As shown at right, UHV pressure increases by about 1.5x from 2 mtorr up to 20 mtorr, an increase of only about 1.8x10⁻¹¹ torr. Therefore, operating the instrument at elevated pressure does not compromise Orbitrap resolution.

5.0x10⁻¹¹ · 4.8x10⁻¹¹ ⋅ 4.6x10⁻¹¹ · 4.4x10⁻¹¹ → 4.2x10⁻¹ 4.0x10 3.8x10⁻¹

Figure 4. The above improvements have positive consequences for the decay of large molecules in the Orbitrap. At right is the decay of the 32+ charge state of carbonic anhydrase, plotted versus IRM pressure. For the "Prototype (Lumos-type)" configuration, decay measurement was not possible above a pressure of 1 mtorr, because only two beats were visible.

In contrast, on the new instrument, decay is nearly constant out to 3 mtorr.





Figure 10. Applying the ΦSDM algorithm specifically to the TMT reporter ions increased the resolution sufficient to baseline resolve TMT isotopologues even when using transients that produce a 30k or 15k resolving power MS2 scan.

(Above) ΦSDM algorithm applied to 0.22 Da windows is sufficient to baseline resolve TMT isotopologues.

(Right) This approach increases the number of identifications for a 50 min gradient with 1 µg of TMT11plex yeast digest standard.

CONCLUSIONS

After modifying hardware and software on a Tribrid mass spectrometer, the instrument's versatility and utility is extended. In particular:

- By raising the IRM pressure and changing the settings for injecting ions into the Orbitrap. sensitivity for higher m/z ions has been increased.
- By integrating the ΦSDM algorithm into the instrument control software, it is now possible to acquire high-resolution spectra in a small percentage of the time required for traditional FFT methods.
- Although TurboTMT is limited to smaller windows in m/z due to available computational hardware, important benefits for TMT experiments are available.

REFERENCES

- 1. Grinfeld et al., Analytical Chemistry 89, 2, 1202-1211 (2016).
- 2. Scheffler et al., ASMS 2016, poster MP312.

ACKNOWLEDGEMENTS

We gratefully acknowledge Mike Goodwin, Chris Mullen, and Raman Mather for helpful discussions.

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.

PO65548-EN0519S

■ 30k TurboTMT ■ 15k TurboTMT

