# Intelligence Driven Metabolomics Workflows: Hardware and Software Innovations for Improved Quantification and Annotation

Susan Bird, Bashar Amer, Rahul Deshpande, Ralf Tautenhahn, Tim Stratton, Brandon Bill, Siegrun Mohring, Sally Webb, and Andreas Hühmer Thermo Fisher Scientific, 355 River Oaks Parkway, San Jose, CA, USA, 95134

#### ORBITRAP-BASED MASS SPECTROMETRY FOR METABOLOMICS

#### High-Quality Data for High-Quality Results

Metabolomics is the comprehensive, qualitative, and quantitative study of all endogenous small molecules in each biological system. Metabolomics is the comprehensive, qualitative, and quantitative study of all endogenous small molecules in each biological system.<sup>1</sup> The collection of these small molecules reflects the biochemical phenotype, which in turn aids in understanding the physiological function and associated pathologies. Yet, the physical-chemical properties of these endogenous compounds are vastly diverse including a range of molecular weight, polarity, structural possibilities, and concentration creating analytical challenges in any metabolimics analysis. The detection of metabolities by electrospray ionization mass spectrometry is further challenged with spectra containing molecular features derived from external sources that are experimentally unrelated or multiple ion species reflecting the same molecule such as adduct formation. <u>Metabolismics analysis requires (1) high resolution to distinguish closely related masses in complex matrices, (2) accurate mass measurements for confident spectral peak assignments. and (3) consistent results from scan-to-scan and run-to-run over extended periods.</u>

Leading Orbitrap-based mass spectrometers (Figure 1) provide high-resolution accurate mass (HRAM) measurements and sensitivity required to measure metabolities in complex matrices. High resolution distinguishes spectral features of similar mass, which is required to differentiate isobaric species and determine line-lisotopic patterns (Figure 2). Accurate mass measurements are paramount for confident spectral assignment (Figure 3).

When combined with advanced separations, high throughput and quantitative capabilities expand the scope of what we know about metabolites and their role in several different areas of study (Figure 1).

Figure 1. Thermo Scientific™ Orbitrap™ based Mass Spectrometers provide HRAM measurements suitable for metabolomics analysis of sample types including animal, plant and cellular components. The Thermo Scientific™ Orbitrap Exploris™ 240 MS instrument is equipped with a quadrupole mass filter and Orbitrap mass analyzer for HRAM MS and MS/MS while the Thermo Scientific™ Orbitrap Tribrid™ Series instruments include a dual pressure linear ion trap in addition to a quadrupole mass filter and Orbitrap mass analyzer for HRAM MS, MS/MS (HCD), CID, UVPD and MS<sup>n</sup> spectra. Optional 1M resolution available a quadrupole mass resolution available.





Thermo Scientific<sup>™</sup> Orbitrap IQ-X Tribrid<sup>™</sup> MS

Figure 2. Fine isotopic distribution for the A1 ion cluster of biotin. Data collection was on an organic extraction of human plasma reference material. NIST SRM 1950, using a Orbitrap Exploris™ 240 MS with a resolution setting of 120K FWHM @ 200 m/z



Creatine C<sub>4</sub>H<sub>9</sub>N<sub>3</sub>O<sub>2</sub> Exact [M+H] m/z 132.0768 122 LINST -0.3 ppm 192 0767 -0.8 ppm 152,0747 182,0767 112-0762 133 2007 122.0767 132.0767

#### Multiple Dissociation for More Spectral Information

Unique to the Tribrid MS platform are multiple fragmentation techniques: (1) higher-energy collision dissociation (HCD) in a high-pressure collision cell (2) collisional-induced dissociation (CID) in an ion trap (3) multi-stage fragmentation or stepwise MS<sup>n</sup> and Ultraviolet Photodissociation (UVPD) for annotating unknown metabolite structures in untargeted metabolomics and lipidomics overvieweite ju 1d (3)

Figure 4. Multiple dissociation techniques provide more fragment ion information for structure characterization and elucidation of unknowns using the Orbitrap IQ-X Tribrid MS. A) CID, HCD, and UVPD dissociation provide unique spec information to distinguish sugar phosphates. B) Higher order fragmentation generates MS<sup>n</sup> spectral trees enabling the systematic breakdown of the flavonoid rutin.



## INTELLIGENCE DRIVEN MASS SPECTROMETRY

AcquireX Data Acquisition to Collect More Biologically Meaningful Data

Insufficient metabolome annotation has limited the biological interpretation of untargeted metabolomics studies. Fragmentation spectra Insufficient metabolome annotation has limited the biological interpretation of untargeted metabolomics sudies. Fragmentiation spectra provide more spectral information to increase the confidence of unknown annotations. The AcquireX intelligent acquisition generates more fragmentation spectra for sample relevant compounds, avoids unrelated background ions, and removes redundancies by using fully automated tetrative inclusion and exclusion lists. Available on the Orbitrap Exploris<sup>112</sup> 240 MS and Tribrid MS systems, AcquireX takes advantage of knowledge to drive acquisition where the blank sample generates a list of ions to exclude for subsequent fragmentation and atrix sample generates a list of frue sample components to prioritize for data dependent MS<sup>2</sup> and MS<sup>5</sup> acquisition (Figure 5). AcquireX leads to information-rich fragmentation of more experimentally relevant compounds

A)

ions 16.21

B)

# MS<sup>2</sup> and MS<sup>n</sup> acquisition (Figure 5). AcquireX leads to inform (Figure 6). Figure 5. The AcquireX Deep Scan acquisition workflow for improved compound annotation.



- First, the AcquireX process obtains the LC-MS data for the blank and a poded sample The AcquireX process creates an exclusion list from the blank and an inclusion list from the sample data. The first data dependent MS<sup>2</sup> run is acquired and the inclusion/exclusion lists are updated after the run On the second injection, MS<sup>2</sup> spectra are acquired for This process is repeated for a user-specified number of intections.

Figure 6. A) Obtaining MS<sup>2</sup> information on compounds vs rygure c. A) Cotaming MS<sup>2</sup> information on compounds vs. background with traditional data dependent acquisition (DDA) compared to AcquireX. B) Sample components with no MS<sup>2</sup> or MS<sup>2</sup> selected for the preferred ion (M + 4) or associated ion (M+ Adduct) for traditional DDA compared to AcquireX, indicating increased MS<sup>2</sup> primarily avoid compound dereplication, during characterization.

AcquireX

beckgroun ions 0.7%

Traditional DDA

#### SOFTWARE BUILT TO ANNOTATE UNKNOWNS

#### Thermo Fisher <sup>™</sup> Compound Discoverer<sup>™</sup> 3.3 Software

Compound Discoverer is a qualitative data-processing application that uses accurate mass data, isotope pattern matching, fragment matching, and mass spectra find the entire product line of Thermo Scientific high-resolution mass spectrometers. The Compound Discoverer 3.3 application includes the following new features:

- New Peak Detection Improvements to mzCloud library search including MS<sup>n</sup> search
- 3. Optimization for large datasets
- Enhancement for GC/MS workflow and molecular networking 4

gure 7. The new peak detection feature can detect aks at very low intensities which would have been





Mass Spectral Library and library match

Mass spectral Library and library is a highly curated, a public library of endogenous and exogenous small molecules containing over <u>5 million fragmentation spectra</u>. Each compound entry in the library generally includes two fragmentation techniques: HCD and CID. This highly diverse spectral library (Table 1) contains true mass spectra generated from purified reference standards. Each spectrum is recalibrated for exact mass and noise removed and is further structurally amotated making this an ultra high-quality spectral library. The mzCloud library is fully integrated into the Compound Discoverer and the Thermo Scientific<sup>™</sup> Mass Frontier<sup>™</sup> spectrum integration spectra endoted from the RXM measurements of the precursor ion, fragmentation spectra privide an additional layer of knowledge about the molecular makeup of a compound and subsequently increase unknown annotation confidence. Fragmentation spectra within the library (Figure 9). Associations based on structural relationships to the parent structure in the reference library enable unknown annotations.





#### SEMI-TARGETED METABOLOMICS

In order to make meaningful biological interpretations of metabolomic data, researchers need two main insights: first, spe identification and quantification of metabolite changes (targeted analysis), and second, an overview of general changes to metabolome (untargeted analysis). Both approaches have pros and cons when using LC-MS, which means researchers s with a choice: a low accuracy overview of total molecular changes, or a detailed yet restricted snapshot of a select group metabolites. archers are often faced t group of

However, a new semi-targeted workflow combining both untargeted and targeted workflows has recently emerged as a middle ground, addressing the limitations of both approaches. The primary focus of the semi-targeted approach is the confident annotation and optionally the accurate quantitation of a targeted set of metabolities, and the secondary focus is to find new molecular connections in a system by performing untargeted analysis <u>on a single injection (i.e.,</u> by reanalyzing (or retro-mining the data).

Figure 10 illustrates the semi-largeted metabolomics workflow utilizing high-resolution accurate mass spectrometry on Orbitrap technology and sophisticated data processing and analysis application solutions for targeted confident (lentification and quantification of metabolites, and untargeted differential analysis and unknown annotation for biomarker discovery. Figure 10, Illustration of the semi-targeted metabolomics workflo



By enabling simultaneous acquisition of hypothesis-led and discovery-led datasets, semi-targeted workflows allow scientists to gain more knowledge about biological systems in a single experiment. Weighing the pros and cons of untargeted and targeted metabolomics has often held back the tremendous potential of metabolomics in life sciences research. Now, the semi-targeted anal looks set to help to unlock this potential.

One of the biggest strengths of semi-targeted metabolomics is the ability to perform targeted and untargeted analysis in a single sample injection. In traditional metabolomics experiments, a sample is injected (analyzed) twice; once for untargeted metabolomics analysis and a second time for targeted analysis. A single injection workflow is particularly advantageous for laboratories that have limited access to samples, time, and resources, and offers a powerful and efficient way to gain more knowledge from valuable biological samples.

#### CONCLUSIONS

· Confidence in compound annotation increases when combining HRAM full scan measurements with multiple dissociations (CID/HCD, MS<sup>2</sup> and MS<sup>4</sup>

Intelligence-driven data acquisition generates more experimentally relevant spectral information and ignores background and redundancies to increase annotation of unknown compounds

Semi-targeted workflow combining both untargeted and targeted workflows to address the limitations of both approaches. The semi-targeted approach ensures the confident annotation and optionally the accurate quantitation of targeted metabolites, in addition, it enables performing untargeted analysis on a single injection.

### REFERENCES

Oliver, S.G., Winson, M.K., Kell, D.B., and Baganz, F. (1998). Systematic functional analysis of the yeast genome. Trends in Biotechnology, 16(9):373-8.

#### TRADEMARKS/LICENSING

© 2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidia This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of



Figure 3. Excellent scan-to-scan mass measurement accuracy was obtained with the Orbitrap Tribrid MS. Subppm mass measurement accuracy for creatine detected over the LC-MS elution profile.

