

Facilitate Biomarker Discovery Using Integrated “Omics” Differential Analysis with High Resolution Accurate LC/MS Approach

Reiko Kiyonami¹, Sergei Snovidia², Devin Drew¹, Rosa Viner¹, Julian Saba¹, David Peake¹, Andreas Huhmer¹, Ken Miller¹,

¹Thermo Fisher Scientific, San Jose, CA, USA; ²Thermo Fisher Scientific, Rockford, IL, USA

ABSTRACT

Application of “Omics” technologies to unbiased clinical biomarker discovery research of various diseases is emerging because comprehensive molecular profiling using these “Omics” technologies could better explain complex disease mechanisms and reach a new level of molecular understanding concerning carcinogenesis, disease progression, and drug-response prediction. Omics approaches are resource-intensive, analytically demanding and require the use of a sophisticated instrument platform and software to identify and quantify hundreds to thousands of molecules with high confidence and accuracy. A complete workflow based on “Omics” profiling also requires well developed sample preparation kits. We have worked on developing multiple “Omics” profiling workflows using high resolution accurate Orbitrap™ mass spectrometers in combination with liquid chromatography separations. In this study, we applied lipidomics, proteomics and glycoproteomics profiling workflows to study of Zucker diabetic fatty (ZDF) and lean rats, a model of type 2 diabetes. Plasma samples collected from three lean and three fatty Zucker rats were analyzed. With the multiple “Omics” workflows developed here we were able to simultaneously identify and quantify hundreds of individual molecular lipid species, proteins and glycopeptides with high confidence and good analytical precision. Significant changes of individual lipid species, proteins and glycopeptides were observed between the two group of rat plasma samples.

INTRODUCTION

HPLC-MS platforms are increasingly used for large scale “Omics” (proteomics, lipidomics, metabolomics, glycomics) experiments to discover potential biomarkers for early disease diagnosis. Each “Omics” data set provides valuable insights into multiple biomarker candidates. However, because properties of a biological system interact with each other, integrating multiple “omics” data will help to understand the system behavior as a whole for unravelling biological regulatory mechanisms to define the emergent properties and help to verify biomarker candidates from each “omics” workflow. Here we conducted lipidomics, proteomics and glycoproteomics studies phenotypic ZDF vs. normal lean rat plasma using a Thermo Scientific™ Orbitrap™ based LC-MS platform. The results from each of the “Omics” workflows are reported.

MATERIALS AND METHODS

Sample Preparation

The rat plasma were purchased from Bioreclamation LLC (Westbury, NY). Rat plasma were recovered from whole blood of Zucker lean (3 lots) and Zucker diabetic fatty (ZDF, 3 lots) using EDTA as anti-coagulant by Bioreclamation. Chloroform, methanol and water were used for the lipid extraction. For proteomics/glycoproteomics experiments, each plasma sample was reduced, alkylated and digested with trypsin, respectively. Tryptic digests (1 mg) from each sample (Lean 1,2,3 and Fat 1,2,3) were labeled with Thermo Scientific™ TMT™ 6-plex (TMTsixplex™) tags, quenched, combined and glycopeptide enrichment was performed using Thermo Scientific™ HyperSep™ Retain AX Cartridges. The flow-through from the cartridge (non-glycopeptides) was collected separately.

Methods

Lipid extracts were analyzed by LC-MS/MS in positive and negative mode on Thermo Scientific™ Dionex™ UltiMate™ 3000 Rapid Separation LC (RSLC) system coupled with a Thermo Scientific™ Q Exactive™ HF hybrid quadrupole-Orbitrap mass spectrometer over a 30 min gradient. Data were collected using a data dependent MS/MS experiment in which a full MS scan (120,000 resolving power FWHM at m/z 200) was followed by top 15 MS/MS scans (30,000 FWHM resolving power at m/z 200). Data analysis was performed using Thermo Scientific™ LipidSearch™ 4.1 software.

Protein digests labeled with TMT sixplex in which the glycopeptides had been depleted were analyzed by nanoflow LC-MS/MS on EASY-nLC™ 1000 coupled with a Thermo Scientific™ Orbitrap Fusion™ Lumos™ mass spectrometer over a 2-hour gradient. Data were collected using top speed data dependent HCD MS/MS experiment in which a full MS scan using 120,000 (FWHM at m/z 200) resolving power was followed by HCD MS/MS scans using 30,000 (FWHM at m/z 200) resolving power with 3 second cycle time. Data analysis were performed using Thermo Scientific™ Proteome Discoverer™ software. Protein annotation was performed using Thermo Scientific™ Protein Center™ software.

Enriched glycopeptides digests labeled with TMT sixplex tags were analyzed by nano flow LC-MS/MS on the Thermo Scientific™ EASY-nLC 1000 coupled with a Thermo Scientific™ Orbitrap Fusion™ Lumos™ Tribrid™ mass spectrometer over a 2-hour gradient. Data were collected using top speed data dependent EThcD MS/MS, followed by SPS MS³ experiments with 3 second cycle time. Mass resolution (FWHM at m/z 200) used was 120,000 for full MS, 30,000 for EThcD MS/MS and 60,000 resolving power for SPS MS³. MS data analysis were performed using Byonic node (Protein Metrics Inc.) in Thermo Scientific™ Proteome Discoverer™ 2.1 software.

RESULTS

Lipidomics

Figure 1. Basepeak Chromatograms of a ZDF Plasma

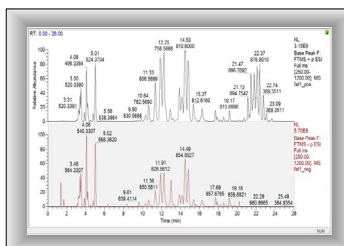
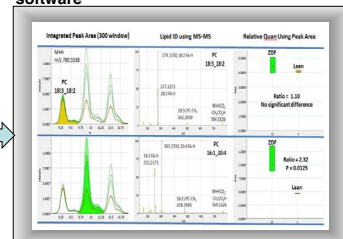


Figure 2. Simultaneous Lipid ID and Relative Quantification Using LipidSearch software



Each plasma sample was analyzed in positive and negative mode, respectively. The C30 reversed-phase HPLC column uniquely offers high shape selectivity and provides improved lipid isomer separation efficiency, yielding higher coverage of the rat plasma lipidome.

All raw files were processed using Lipid Search™ 4.1 SP1 software. The ultra high resolution offered by the Q Exactive HF MS enables very narrow mass search windows (3 ppm for MS and 5 ppm for MS/MS) being used for database searching, providing high confidence lipid identification and precise quantification results.

Table 1. Summary of Lipid IDs and Quantitation from the Rat Plasma

Class	Filtered*	p < 0.05
PA	1	0
PC	135	93
LPC	110	77
PE	39	0
LPE	4	3
PG	4	2
PI	42	30
PS	3	3
Car	21	13
CerG1	13	10
SM	28	16
AcCa	19	14
ChE	11	5
StE	5	3
StE	3	0
DG	30	29
TG	2083	522
Total	750	349

The lipid annotations for each individual sample are aligned within a time window and the results are merged into a single report. Positive and negative ion adduct annotations are grouped for lipids eluting at the same retention time. Accurate-mass extracted chromatograms are integrated for each identified lipid precursor ion and peak areas are obtained for quantitation. t-Tests and calculated p-values determine which molecular species are significantly different between sample and control groups. 750 Molecular lipid species across 17 lipid classes were identified after filtering. Fold-changes of the identified lipid species between the lean plasma and the diabetic plasma groups were determined. Significantly different concentration changes (p < 0.05) were observed for 349 lipid species.

The figure consists of three vertically stacked bar charts, each comparing two cases: S1: ZDF (green bars) and C: Lean (orange bars). The y-axis for all charts represents a value, with scales varying between charts. The x-axis for each chart has two categories, labeled with symbols that appear to be σ and π .

- Top Chart:** The y-axis ranges from 0.00E+00 to 2.50E+11. For the σ category, the C: Lean bar is approximately 2.00E+11 and the S1: ZDF bar is approximately 2.50E+11. For the π category, the C: Lean bar is approximately 2.00E+11 and the S1: ZDF bar is approximately 2.50E+11.
- Middle Chart:** The y-axis ranges from 0.00E+00 to 3.50E+11. For the σ category, the C: Lean bar is approximately 2.50E+11 and the S1: ZDF bar is approximately 3.00E+11. For the π category, the C: Lean bar is approximately 2.50E+11 and the S1: ZDF bar is approximately 3.00E+11.
- Bottom Chart:** The y-axis ranges from 0.00E+00 to 7.00E+09. For the σ category, the C: Lean bar is approximately 2.00E+09 and the S1: ZDF bar is approximately 1.00E+09. For the π category, the C: Lean bar is approximately 2.50E+09 and the S1: ZDF bar is approximately 2.00E+09.

Lipidomics

Proteomics/Glycoproteomics

Wiki Pathways

Wilda Pathways table

Adipogenesis
Blood Clotting Cascade
Brain derived neurotrophic
Complement Activation, Complement and Coagulation
Fetal Adhesion
Inflammatory Response
Myometrial Relaxation and Con...
Tet...

Unannotated (28)
Steroid Biosynthesis (WP66) (1)
Regulation of Actin Cytoskeleton (12)

high zdf

General Data Information
Source: Ensembl (14/04/2012)

high zdf
Reference data set

Over-represented Wiki pathways

Analysis data set: high zdf Reference data set: Rattus norvegicus (G1)

Description	Occurrence	Count	Ref. Count	Raw p-value	FDR p-value
Complement and Coagulation Cascades (WP347)		14	274	1.68E-18	1.51E-18
Blood Clotting Cascade (WP255)		7	78	2.57E-11	1.16E-10
Complement Activation, Classical Pathway (WP1)		3	77	2.68E-4	8.04E-4
Inflammatory Response Pathway (WP40)		2	127	1.87E-2	4.21E-2

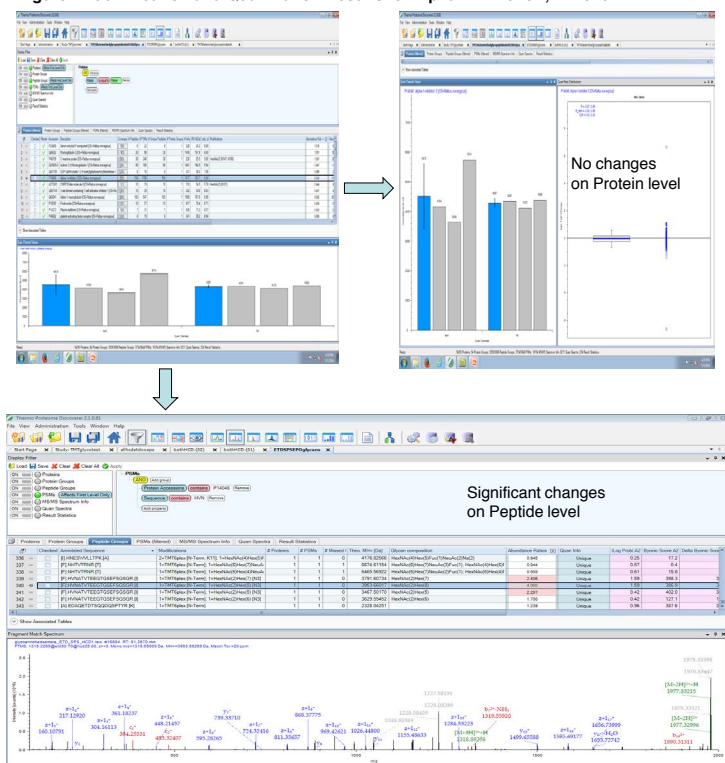
#191 61 members - sum=70.16

KEEVAWTELLAKI
KESTLEPSEDCLYLNYSPAD...
RLGMLPSTIGERHG
RLSPREVENTIONALFINK...
PPPPVQLVLTYSNEN
RLVDEKQDPINQVFXA
KYVLESGTQPVAVLGPFPK...
KINKEOCQDLKADORA
KPLEVNNYHSEAFVNFADSE...
[52 more]

AQR7 AQR8 AQR9 AQR1 AQR2 AQR3 AQR4 AQR5 AQR6

[illegible]

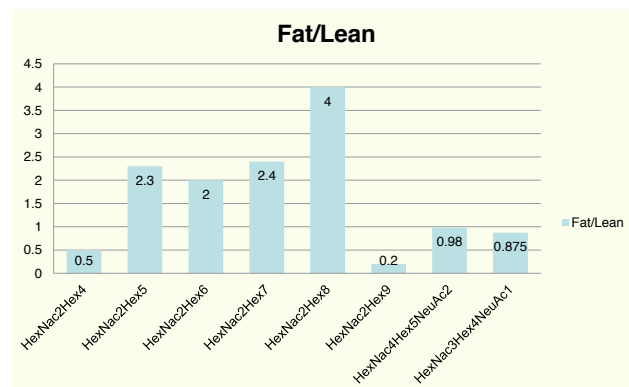
Figure 7. Identification and Quantitation Results for Alpha-1 inhibitor , P14046



The protein of the Alpha-1 inhibitor did not show changes in protein levels. However, significant changes in glycopeptide profiles were observed as shown in Figure 7. For example, the amount of high mannose glycoforms for site N321 increased 2-4 times in ZDF rat plasma (Figure 8), but there are no changes for hybrid/complex glycosylated species.

Proteomics/Glycoproteomics

Figure 8. Expression Fold Changes of Peptide HVNATVTEEGTGSEFGSGR- Site N321 in Alpha-1 inhibitor, P14046



The up and down changes in the peptide level with different glycan composition. High mannose species are up and complex/hybrid glycans are down.

CONCLUSIONS

- Complete multiple "Omics" workflows were developed on high resolution accurate mass Orbitrap – based mass spectrometers coupled with liquid chromatography separation and sophisticated data processing software. The ultra high resolution offered by Orbitrap detector enables another dimension of mass separation, yielding confident unknown molecular identification and precise quantification.
- 750 Molecular lipid species were identified and quantified from the rat plasma samples. The relative concentration changes between the ZDF and lean plasma groups were measured at both total lipid class composition and individual lipid molecular species composition, providing more potential biological insights into understanding the complex processes that lead to the development of diabetes.
- Unique data dependent ETHcD MS/MS and SPS MS³ workflows enabled very precise protein/peptide quantitation in a high throughput fashion. Unambiguous glycopeptide characterization was performed using HRAM ETHcD spectra. Precise relative quantitation was carried out using HRAM SPS HCD MS³ data.
- Significant increases of glycopeptides containing high mannose species were observed between the ZDF and lean plasma without any changes at the protein level. Further studies will be done to see if these glycan composition changes are linked with lipid molecular species changes.

www.thermofisher.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. Lipid Search is a registered trademark of MKI. TMT and TMTsixplex are trademarks of Proteome Sciences plc. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others. 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 +43 1 333 50 34 0
Australia +61 3 9757 4300
Austria +43 810 282 206
Belgium +32 53 73 42 41
Brazil +55 11 2730 3006
Canada +1 800 530 8447
China 800 810 5118 (free call domestic)
 400 650 5118

Denmark +45 70 23 62 60
Europe-Other +43 1 333 50 34 0
Finland +358 10 3292 200
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9494
Italy +39 02 950 591

Japan +81 6 6885 1213
Korea +82 2 3420 8600
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
Norway +46 8 556 468 00

Russia/CIS +43 1 333 50 34 0
Singapore +65 6289 1190
Sweden +46 8 556 468 00
Switzerland +41 61 716 77 00
Taiwan +886 2 8751 6655
UK/Ireland +44 1442 233555
USA +1 800 532 4752
 PN64735-EN 0616S

Thermo
 SCIENTIFIC
 A Thermo Fisher Scientific Brand