

Processing of a Complex Lipid Dataset for the NIST Inter-laboratory Comparison Exercise for Lipidomics Measurements in Human Serum and Plasma

David A. Peake¹, Reiko Kiyonami¹, Yasuto Yokoi², Yukihiro Fukamachi², Ken Miller¹

¹Thermo Fisher Scientific, San Jose, CA, USA; ²Mitsui Knowledge Industry, Tokyo, Japan

Introduction

Lipids play a key role in cell, tissue and organ physiology with diseases such as cancer and diabetes which involve disruption of their metabolic enzymes and pathways. Identification of unique lipid biomarkers to distinguish healthy humans compared to those with a disease can have an impact on the early detection of diseases and personalized medicine.

Identification of lipids by untargeted lipidomics requires sophisticated software with an extensive lipid database. In addition, the mass spectrometer employed must be capable of separating many overlapping isobaric and isomeric lipid ions. We present here the details and challenges of the data processing of NIST plasma and serum extracts using the latest version of LipidSearch software. New algorithms were introduced specifically to reduce false positives and to automate the data review.

Thermo Scientific™ LipidSearch™ software was used for lipid identification through a database search of precursor accurate masses and their predicted fragment ions. Each lipid identification is ranked by mass tolerance, match to the theoretical fragment ions and fraction of total MS² intensity. The number of lipid species identified in each LC-dd-MS² experiment were assessed at sum composition (MS) and isomer (MS²) levels. Potential lipid species were identified using the predicted MS-MS fragments for molecular species observed in positive or negative ion mode.

The data for each run were aligned within a chromatographic time window and positive and negative ion annotations were merged into the results table. This approach provides lipid annotation that reflects the appropriate level of MS² fragment ions from the complete dataset giving higher confidence in lipid identifications. The merged results were filtered by main adduct ion, ID quality, signal-to-noise, peak area and relative standard deviation; manual integration was performed if necessary prior to estimating concentration relative to an internal standard for each lipid class. These results demonstrate that in a 60 min LC-MS run that it is possible to identify and quantify approximately 1000 isomeric lipid species from human plasma using an experimental C30 UHPLC column.

Methods

LC-MS Sample Preparation. Aliquots of human plasma or serum (80µL, Table 1) were extracted using the method of Bligh and Dyer.

LC-MS and dd-MS² Method. A Thermo Scientific™ Dionex™ UltiMate™ 3000 Rapid Separation LC (RSLC) system and Thermo Scientific™ Q Exactive HF™ instrument were employed as described (1). The HPLC separation was achieved with a 2.1 x 250mm, 1.9µm, C30 prototype column and the MS analysis was performed at 120K resolution and MS² at 30K resolution (FWHM at m/z 200).

Data Analysis. The data were processed using a HP Z840 workstation equipped with dual Hex-core Xeon processors (E5-2643v3, 2133 6C, 3.4 GHz), 64 GB of RAM (DDR4-2133), and a 500GB solid state hard drive.

Data Processing

LipidSearch software was used for lipid identification and relative quantitation using the workflow shown in Figure 1 (2, 3). LC-MS data (24 raw files, 5.9 GB total) containing MS and data dependent-MS² were searched using the parameters listed in Table 2. For each MS² spectrum, search results are summarized for lipid species matching the predicted fragmentation pattern from the database with match score and occupancy indicating the fit (Figure 2). The average number of sum composition lipids identified from LC-MS runs collected in positive ion or negative ion mode was 211 and 867, respectively. The same lipid annotations within ± 0.1 min were merged into the aligned results (Figure 3). The total number of lipids identified in the entire data set are summarized in the Table 3. Lower quality annotations and false positives were removed using a combined set of data filters prior to estimation of concentration relative to an internal standard for each class.

TABLE 1. NIST Human Plasma/Serum Samples

| Group | NIST ID | Sample Type | Sample Description |
|-------|----------|--------------|--|
| C | SRM 1950 | Human plasma | Equal number of men and women |
| S-1 | 2378-1 | Human serum | Donors took fish oil supplements |
| S-2 | 2378-2 | Human serum | Donors took flaxseed oil supplements |
| S-3 | 2378-3 | Human serum | Donors did not take fish or flaxseed oil |

TABLE 2. LipidSearch Conditions

| Search Parameter | Settings | Units |
|---------------------------|--|-------|
| Precursor mass tolerance | 5.0 | ppm |
| Product mass tolerance | 5.0 | ppm |
| Prod. Intensity threshold | 1.0 | % |
| m-Score threshold/display | 2.0 / 5.0 | |
| Quan m/z tolerance | +/- 5.0 | ppm |
| Quan range | +/- 1.0 | min |
| Main isomer peak filter | ON | |
| ID Quality filter | A,B,C,D | |
| Adducts (pos ion) | H, NH ₄ , Na | |
| Adducts (neg ion) | -H, -2H, +CH ₃ CO ₂ | |
| Lipid Sub-Classes | Phospholipids: LPC, PC, LPE, PE, LPS, PS, LPG, PG, LPI, PI, LPA, PA, CL Sphingolipids: SM, Cer, CerG1, CerG2, CerG3 Glycerolipids: DG, TG Neutral lipids: ChE, Co | |
| Merge Parameter | Settings | Units |
| Retention time tolerance | 0.20 | min |
| All isomer peak filter | ON | |
| m-Score threshold | 5.0 | |
| ID Quality filter | A,B,C, D | |

Results

LC-MS/MS Data Processing Workflow using LipidSearch software (Figure 1).

- 1) Peak Detection.** Read raw files, MSⁿ and precursor ion accurate masses.
- 2) Identification.** Candidate molecular species are identified by searching a large database > 10E+7 entries of accurate masses (lipid precursor and fragment ions) predicted from each potential lipid structure and positive/negative ion adducts.
- 3) Alignment.** The search results for each individual sample are aligned within a time window and the results are merged into a single report.
- 4) Quantification.** The accurate-mass extracted ion chromatograms are integrated for each identified lipid precursor and the peak areas are obtained. Analyte concentration for each lipid class was calculated relative to the concentration of internal standard.
- 5) Statistical Analysis.** t-Tests determine which lipid species are significantly different between sample vs. control groups, and results are displayed in a whisker plot.

FIGURE 1. LipidSearch Workflow.

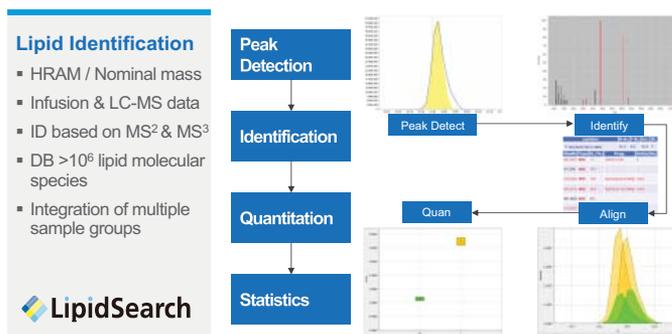


FIGURE 2. LipidSearch Results for Neg. Ion m/z 762.5289, Rt = 18.05 min.

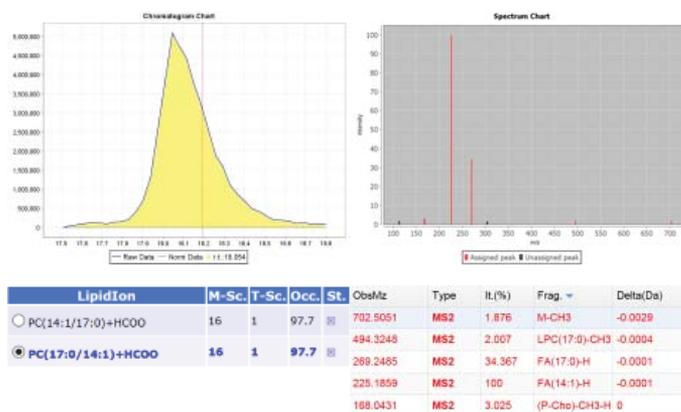


FIGURE 3. Aligned Peak Areas for PC Internal Standard 17:0/14:1, M+H⁺.

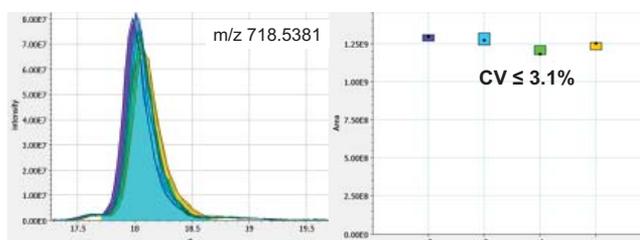


TABLE 3. Number of Lipid Species Identified after Merging Data, Filtered IDs, and Number of Lipid Species Reported and Quantified for Each Lipid Sub-class.

| Lipid Class | No. Species Unfiltered | No. Species Filtered | No. Species Reported |
|--------------|------------------------|----------------------|----------------------|
| ChE | 22 | 19 | 18 |
| DG | 81 | 46 | 45 |
| TG | 665 | 468 | 452 |
| PC | 508 | 221 | 220 |
| LPC | 106 | 58 | 57 |
| PE/LPE | 53 | 33 | 32 |
| PI | 34 | 25 | 24 |
| Cer | 33 | 22 | 21 |
| CerG | 20 | 13 | 13 |
| SM | 138 | 92 | 91 |
| Total | 1660 | 997 | 973 |

Figure 4 shows the relative levels of four different TG 54:6 isomers in the human serum and plasma samples: 1) TG(18:1/18:2/18:3) at 40.04 min, 2) TG(18:2)₃ at 40.25 min, 3) TG(16:0/18:1/20:5) at 41.22 min and 4) TG(16:0)₂(22:6) at 41.90 min, retention time. Serum 2 (donors taking flaxseed oil supplement) is significantly higher in triglyceride containing 18:2 and 18:3 (ALA) fatty acids whereas Serum 1 (donors taking fish oil supplement) contains more TG with polyunsaturated 20:5 (DPA) and 22:6 (DHA).

FIGURE 4. Comparison of Isomeric TG 54:6 Species in Human Serum/Plasma.

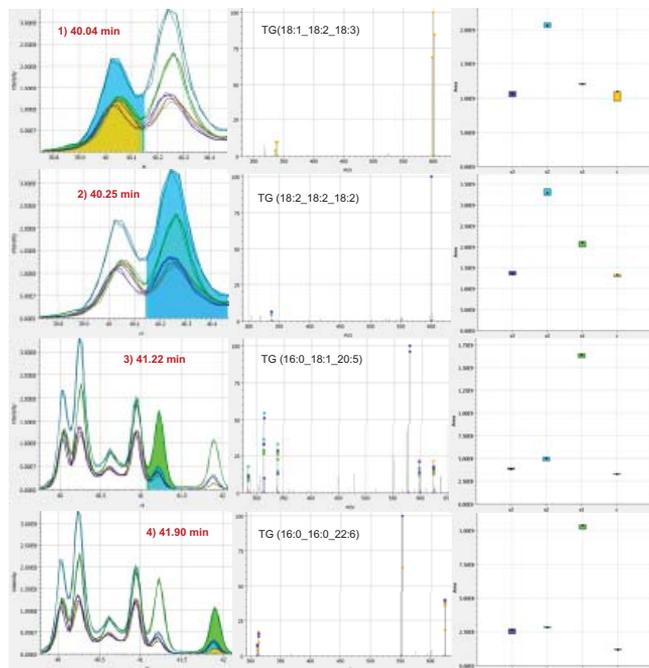


TABLE 4. Concentration of TG 54:6 in NIST Human Serum and Plasma Samples.

| TG Species | Rt, min | Conc., nmol/mL | | | | CV, % | | | |
|---------------------|---------|----------------|------|------|------|-------|-----|-----|-----|
| | | S-1 | S-2 | S-3 | C | S-1 | S-2 | S-3 | C |
| TG (18:1/18:2/18:3) | 40.04 | 3.22 | 5.31 | 2.26 | 3.01 | 3.1 | 2.6 | 2.4 | 5.1 |
| TG (18:2/18:2/18:2) | 40.25 | 5.57 | 8.49 | 2.91 | 3.79 | 4.1 | 3.5 | 1.7 | 3.0 |
| TG (54:6) | 40.62 | 3.06 | 3.31 | 1.50 | 2.10 | 3.2 | 5.5 | 7.3 | 6.0 |
| TG (54:6) | 40.94 | 4.79 | 4.84 | 2.43 | 3.37 | 4.4 | 4.1 | 6.6 | 4.1 |
| TG (16:0/18:1/20:5) | 41.21 | 4.38 | 1.28 | 0.82 | 0.95 | 1.6 | 5.1 | 2.7 | 4.7 |
| TG (16:0/16:0/22:6) | 41.90 | 2.76 | 0.73 | 0.55 | 0.34 | 2.2 | 2.1 | 6.1 | 1.1 |

Conclusion

- LipidSearch software provides an automated workflow for processing high quality Orbitrap LC-MS/MS untargeted lipidomics data and enables reliable and comprehensive lipid identification and quantification.
- The high mass accuracy in both MS (120K) and MS² (30K) obtained with the Q Exactive HF instrument allows confident lipid species identification from the highly complex human serum and plasma extracts. Almost a thousand of species are identified automatically and quantified from a single LC-MS run.
- Orbitrap data combined with LipidSearch software allows simultaneous lipid ID with high coverage and quantitation. Each lipid ID was obtained with a single high quality Orbitrap MS² scan over 4 orders of concentration dynamic range.
- The CV of the technical replicates for a majority of the quantified lipids were typically below 15% showing the excellent reproducibility of the Q Exactive HF.

References

1. R Kiyonami, D A Peake, X Liu and Y Huang "Large Scale Lipid Profiling of a Human Serum Lipidome Using a High Resolution Accurate Mass LC/MS/MS Approach" presented at the LIPID MAPS Annual Meeting 2015, May 12-13, 2015, La Jolla, CA.
2. R Taguchi, *et al.*, *J. Chrom. A*, **2010**, *1217*, 4229–4239. doi: 10.1016/j.chroma.2010.04.034.
3. T Yamada, *et al.*, *J. Chrom. A*, **2013**, *1292*, 211-218. doi: 10.1016/j.chroma.2013.01.078.

www.thermoscientific.com

©2015 Thermo Fisher Scientific Inc. All rights reserved. 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 manners 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
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 45 453 9100
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
Spain +34 914 845 965
Sweden +46 8 556 468 00
Switzerland +41 61 716 77 00
UK +44 1442 233555
USA +1 800 532 4752

Thermo
 SCIENTIFIC

A Thermo Fisher Scientific Brand