

Software Utilizing Positive and Negative Ion MS²/MS³ HCD and CID Spectra for Improved MSⁿ Lipid Identification

David A Peake¹, Reiko Kiyonami¹, Daniel Gachotte², Gavin E Reid³, Yasuto Yokoi⁴, and Andreas Hühmer¹

¹Thermo Fisher Scientific, San Jose, CA, United States; ²CORTEVA Agriscience, Indianapolis, IN, United States; ³The University of Melbourne, Parkville, Victoria, Australia; ⁴Mitsui Knowledge Industry, Atago, Tokyo, Minato-ku, Japan

ABSTRACT

Thermo Scientific™ LipidSearch™ software version 4.2 for untargeted lipidomics with LC dd-MSⁿ analysis provides enhanced lipid characterization by combining information obtained in HCD and CID MS² and MS³ product ion spectra. This lipid workflow is demonstrated on a new Tribrid™ Orbitrap™ mass spectrometer designed for characterization of small molecules such as metabolites and lipids. Total lipid extracts from western corn rootworm larvae were analyzed giving high confidence in lipid annotations and good reproducibility for relative and estimated quantitation of almost one thousand lipid species. Phosphatidylcholine and triglyceride lipids were automatically targeted by specific product ion neutral losses, which enabled further CID MS²/MS³ characterization during the same dd-MSⁿ acquisition cycle. Functional lipids such as PI isomers and sterol esters are readily annotated and quantified at low levels. Different phenotypes are observed between two isomeric fatty acyl combinations: PI 18:1_18:3 and PI 18:2/18:2.

INTRODUCTION

Lipidomics is a rapidly growing field of study that is crucial for understanding the underlying mechanisms for disease progression. The application of untargeted lipidomics profiling to phenotypic analysis of a wide range of plant tissues and organisms is becoming a more important aspect of agricultural research. Identification of lipids by untargeted lipidomics requires sophisticated software with an extensive lipid database. We present here the data processing of western corn rootworm (WCR) lipid extracts using the latest version of LipidSearch software. New algorithms were introduced specifically to reduce false positives, improve quantitation using labeled internal standards, and automate searching of MSⁿ data obtained by higher-energy collisional dissociation (HCD) and linear ion trap collisional induced dissociation (CID) fragmentation methods.

MATERIALS AND METHODS

Total lipid extracts from western corn rootworm larvae¹ were separated using a Thermo Scientific™ Accucore™ 2.1 x 150 mm C30, 2.7 μm column and Thermo Scientific™ Vanquish™ chromatograph. LC-MSⁿ analysis was performed using a Thermo Scientific™ Orbitrap ID-X™ Tribrid™ mass spectrometer (Figure 1) using Acquire-X™ experimental workflow for lipid characterization (Figure 2). Lipid extracts were spiked with SPLASH[®] LipidoMIX[®] (Avanti Polar Lipids, Inc.) internal standards to estimate the concentration of annotated lipid species.

LC/MS at 120K resolution (FWHM @ m/z 200) and data-dependent HCD MS² experiments (15K resolution) were performed in positive and negative ion modes. During each 1.5 s cycle of the untargeted dd-MSⁿ profiling method, additional targeted product ion (m/z 184.0733) or neutral loss (fatty acid + NH₃) CID MS² and MS³ experiments were selectively triggered to provide higher quality characterization of phosphatidylcholine (PC) and triglyceride (TG) lipids. Prior to analysis of the biological samples, LC/MS (120K resolution) analyses of blank and pooled samples were used to automatically create an exclusion list for background ions and inclusion list for sample-relevant ions. Both lists were dynamically updated following subsequent LC/MSⁿ analyses of biological replicates.

LC/MSⁿ datasets were processed using LipidSearch 4.2.9 beta software (Figure 3) with an expanded lipid database (Table 1), improved peak detection during alignment of annotated peaks, and better rejection of false positives (Table 2). Multiple HCD MS², CID MS², and MS³ product ion mass spectra for the same precursor ions were automatically combined to provide more comprehensive annotation for hundreds of lipid molecular species. The data processing parameters used for the WCR lipid samples (Figure 4) are shown in Table 3.

RESULTS

LipidSearch software was used to search MS² and MS³ spectra against the predicted product ions and neutral losses for all potential lipid species within precursor and product ion mass tolerances. For each LC-dd-MSⁿ analysis, potential lipid species were identified separately from positive or negative ion adducts. The data for each biological replicate were aligned within a chromatographic time window by combining the positive and negative ion annotations and merging these into a single lipid annotation in the results table. This approach provides lipid annotations that reflect the appropriate level of MS²/MS³ product ions and neutral losses from the entire dataset giving higher confidence in lipid identifications.

The alignment results were filtered by minimum number of data points, signal-to-noise ratio, intensity ratio, main adduct ion, and ID quality. Compared to results generated only from dd-MS² HCD data, the combination of HCD and CID MS²/MS³ gives significantly higher quality lipid identifications for PC and TG lipids in the same analysis time. We will continue to refine this approach based on published lipid MS/MS compilations.²

LipidSearch software version 4.2 has been modified to use the shorthand notation suggested by Liebisch, *et al.*³ and sphingolipid annotations now reflect the number of hydroxyl groups (m = monohydroxy, d = dihydroxy, and t = trihydroxy), the number of carbons in the long-chain base, and the number of double bonds. For example, d18:1/18:0 ceramide becomes Cer[d18:1_18:0] and sphinganine as SPH[d18:0]. To avoid confusion regarding very long chain ceramides found in stratum corneum lipid extracts with a very different nomenclature,⁴ these lipids are now searched using a separate skin lipids target database.

The concentration of lipid species with a deuterated standard were estimated using the internal standard concentration times the peak area ratio of analyte to internal standard (Table 4). The number of lipid species per sub-class after filtering rejected peaks (annotation grades A+B) are summarized in Table 5.

Figure 1. Untargeted lipidomics high-resolution LC-MSⁿ workflow using C30 reversed-phase separation, Orbitrap ID-X Tribrid mass spectrometer, and LipidSearch software

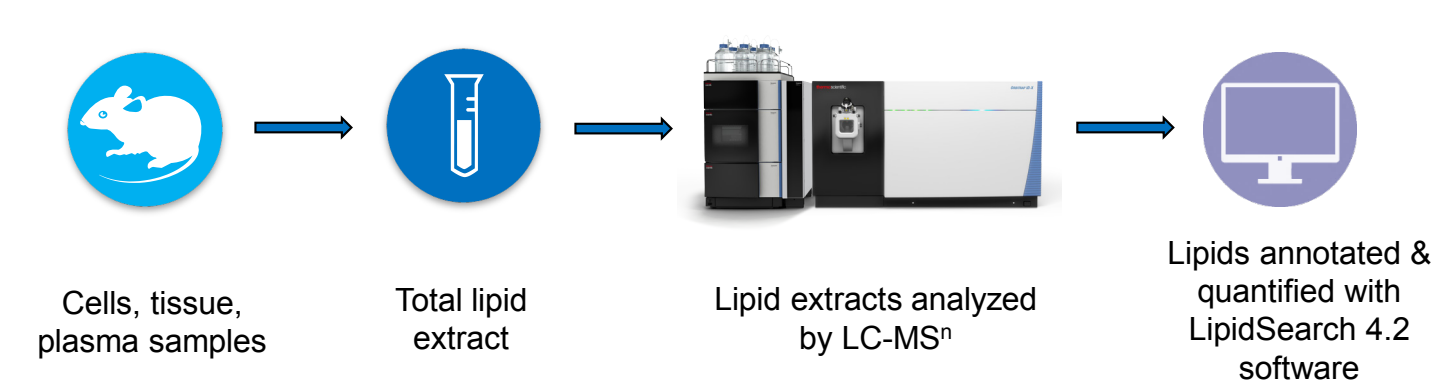


Figure 2. Untargeted LC-MSⁿ (with/without inclusion) followed by targeted CID MS² and MS³

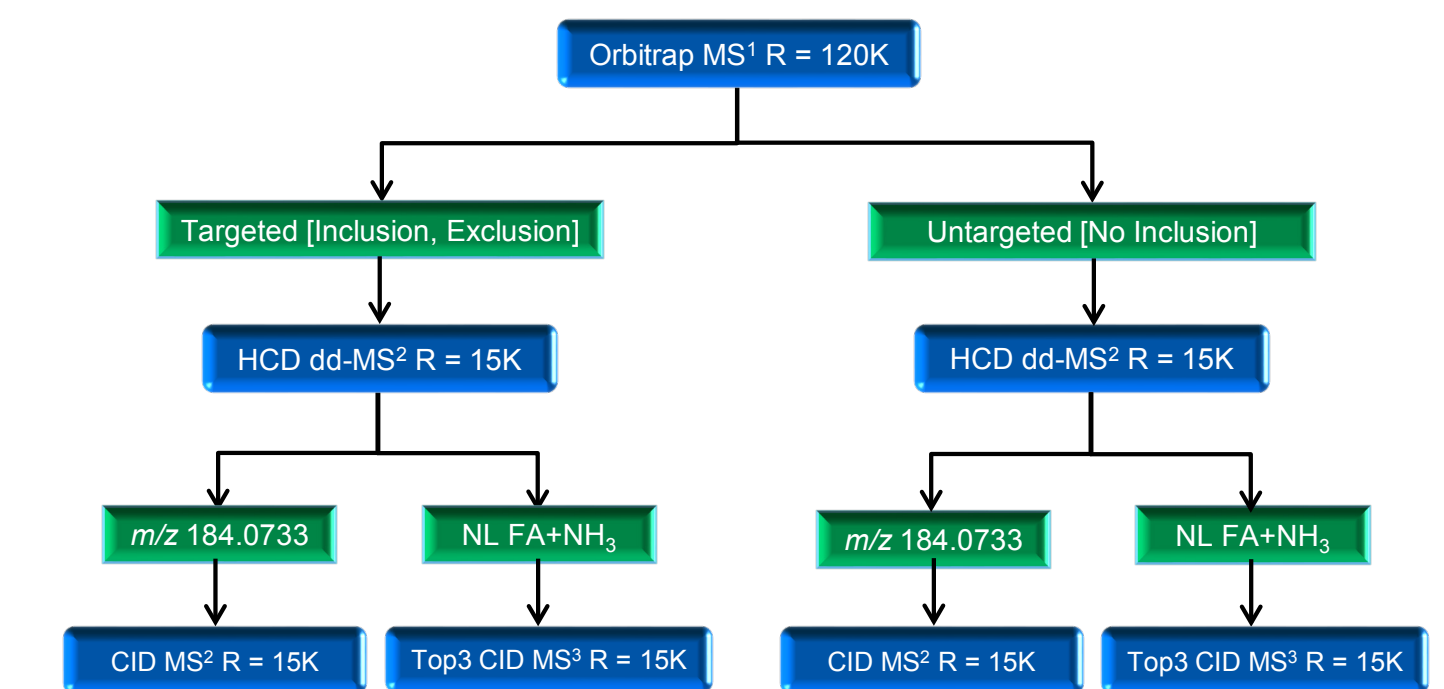


Figure 3. LipidSearch software for lipid annotation and quantitation for LC-MSⁿ data

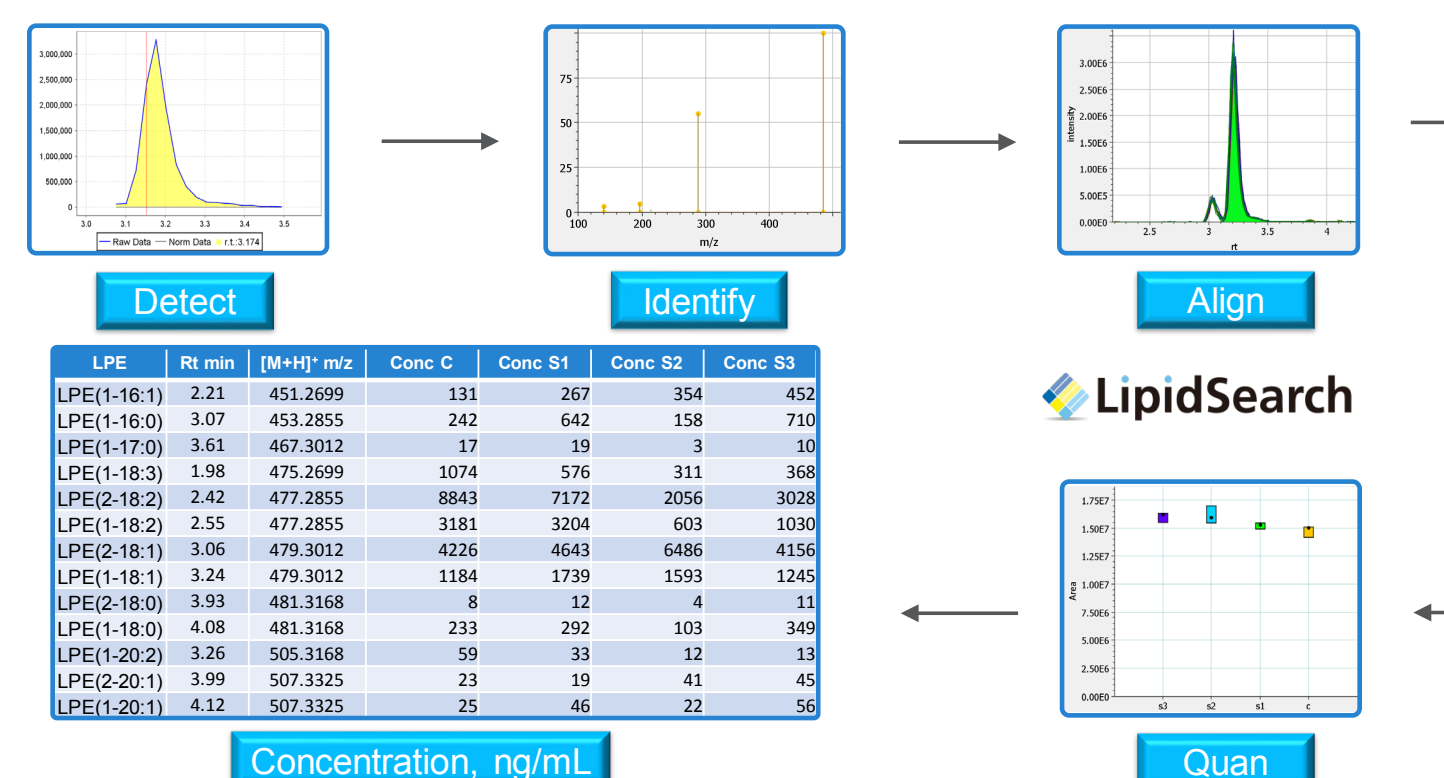


Figure 4. Western corn rootworm larvae lipid extract at two different growth stages and diets

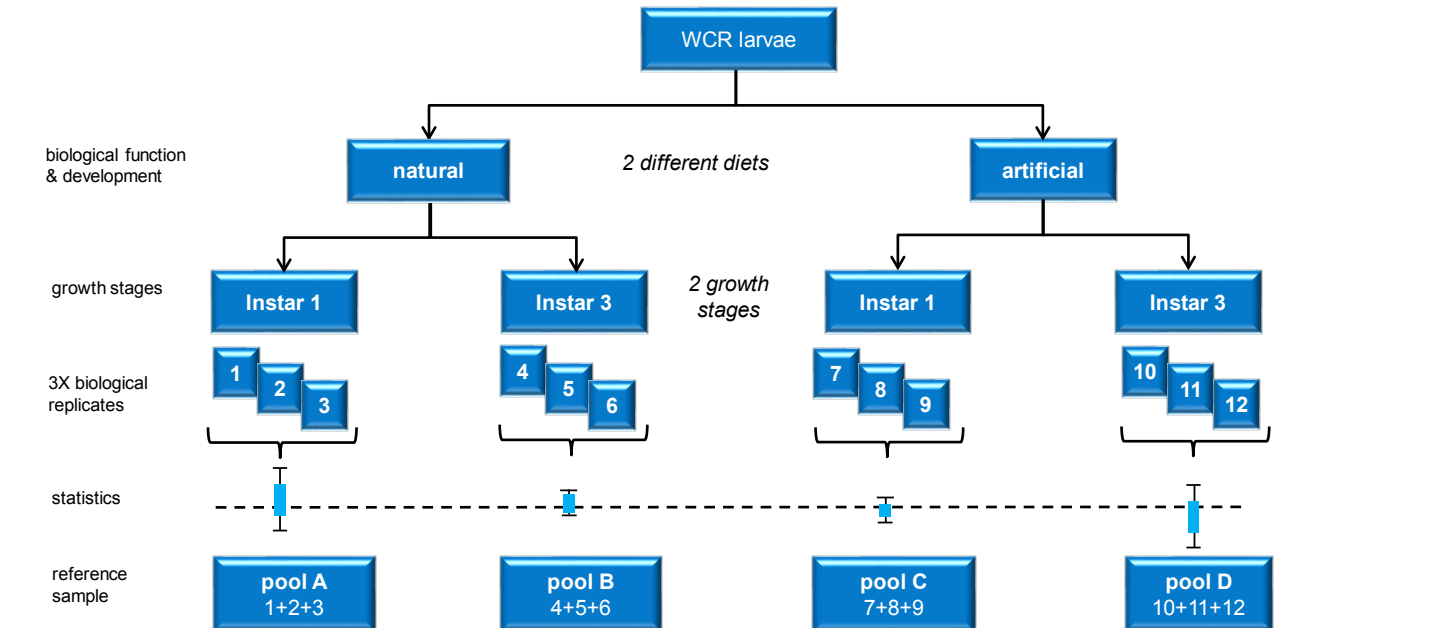
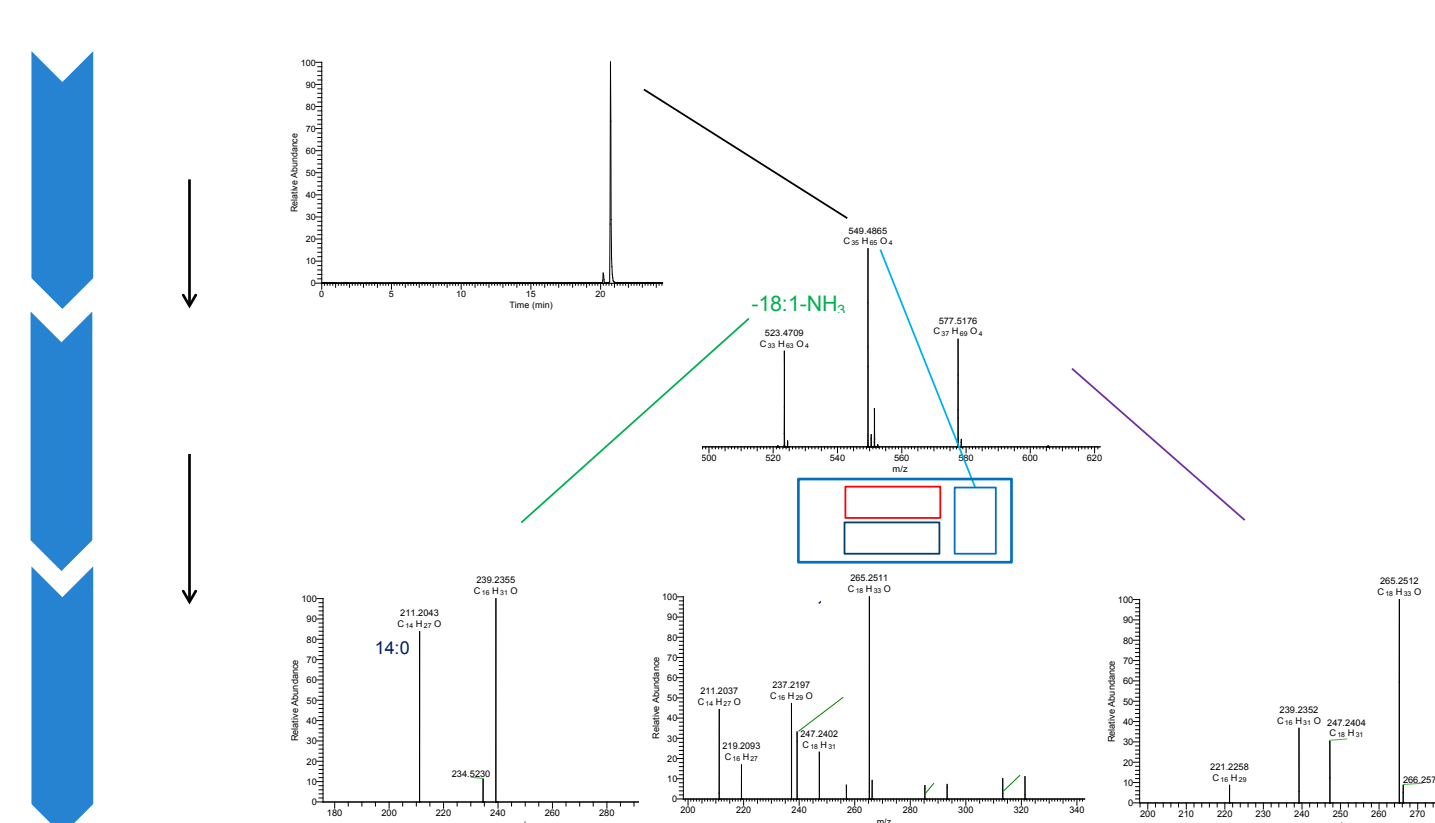


Figure 5. HCD MS² and CID MS³ spectra of Triacylglycerol 48:1 [M+NH₄]⁺ at 20.66 min



Triglyceride ammonium adducts dissociate in MS² to give up to 3 product ions formed by the loss of fatty acids. MS³ of these product ions may correspond to several different fatty acyl combinations. In Figure 5, TG 48:1 gives major product ions corresponding to loss of 18:1 (green), 16:0 (blue), and 14:0 (purple) fatty acids. The MS³ spectra of 882.7541 → 523.4709 (NL 18:1+NH₃) forms acyl product ions for 14:0 and 16:0, giving a 14:0_18:1_16:0 annotation. Similarly, 882.7541 → 577.5176 NL of 14:0+NH₃ produces 16:0, 18:1 giving the same 14:0_18:1_16:0 annotation. However, 882.7541 → 549.4865 NL of 16:0+NH₃ gives a mixture of product ions: 14:0, 18:1 and 16:0, 16:1 fatty acyl ions. Thus, the TG 48:1 is a mixture of TG 14:0_18:1_16:0 and 16:0_16:1_16:0 isomers. This example demonstrates that MS³ enables confident annotation of co-eluting triglyceride isomers typically found in complex total lipid extracts of biological origin.

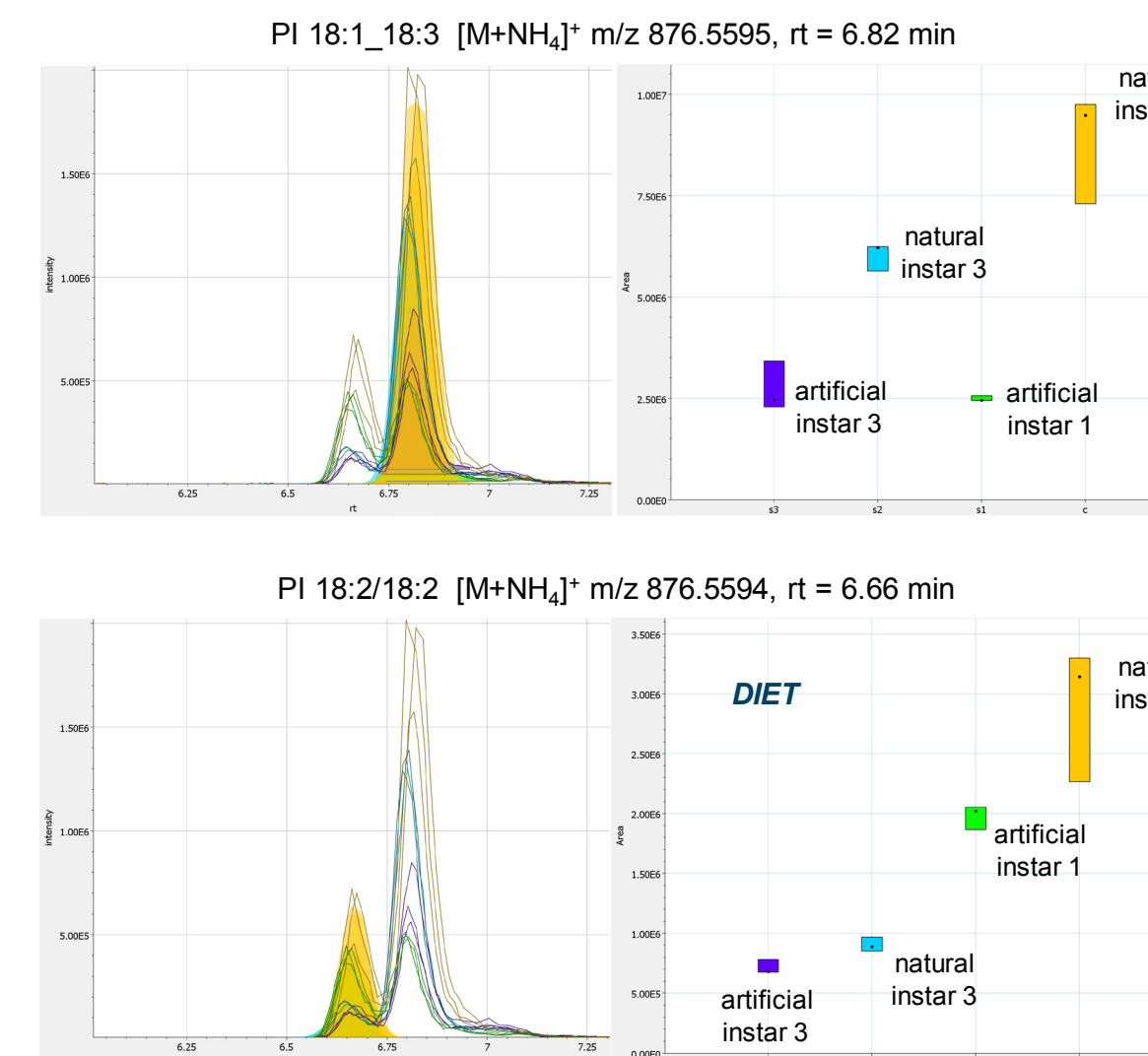
Table 1. Database and lipid nomenclature modifications in LipidSearch 4.2 software

Modification	Description
Added AcHexSIE, AcHexSIE, AcHexZyE, AcHexCmE, AcHexCHE	Acylhexosyl sterol esters
Added bis-methyl LPA, PA, PC, PE, PG, PS	Methylation of phosphate for quantitation
Added methyl PC	Methylation of phosphate for quantitation
Added MLCL, DLCL	Mono-lyso and di-lyso cardiolipins
Added SL	Seminalipids (glycerol analog of sulfatide)
Added 15:0-18:1(d) DG, PA, PC, PE, PG, PI, PS; Chol(d)	SPLASH internal standards for quantitation (Avanti Polar Lipids, Inc.)
Added CerPE	Ceramide phosphatidylethanolamine
Added NAE	N-acyl ethanolamine
Added WE	Wax esters of fatty acids
Modified sphingolipid nomenclature	As proposed by G. Liebisch, <i>et al.</i> ³
Created skin ceramide database	Very long chain skin ceramides ⁴
Added FA chloride adducts	Negative ion APCI for stratum corneum lipids ⁴
Added [M+H-H ₂ O] ⁺ and [M-CH ₃] ⁺ adducts	Detection of cholesterol and phosphatidylcholine in-source product ions

Table 2. New features and improvements in LipidSearch 4.2 software

Feature/Improvement	Description
	Minimum number of data points
	Peak intensity ratio threshold
	Improved signal-to-noise ratio threshold
	Filter duplicate IDs at the same retention time
	Automatic filtering of rejected peaks
	Combines HCD MS ² with targeted (NL or product ion) CID MS ² /MS ³ spectra for enhancing annotation of PC, TG, and other lipid species
	MS ⁿ Num – reports the number of annotations that include CID MS ³
Improved confidence in annotation by combining HCD and CID MS ⁿ searches	Quantification expert using SPLASH internal standard concentration
Estimated quantitation added	Quantification expert using SPLASH internal standard concentration
Broader lipid sub-class coverage	Expanded lipid database adds 26 more lipid sub-classes (total 92)

Figure 6. Low-abundant PI isomers show phenotypical differences



Low-abundant lipids (e.g., PI 36:4) gave significant differences between growth stages: PI 18:1_18:3 (m/z 876.5595, 6.82 min) and between dietary conditions: PI 18:2/18:2 (m/z 876.5594, 6.66 min).

Table 3. LipidSearch 4.2.9 beta software parameters

Search Parameter	Setting	Units
Precursor mass tolerance	5	ppm
Product mass tolerance	10	ppm
Prod. intensity threshold	1.0	%
m-Score threshold/display	2.0/5.0	
Quan m/z tolerance	+/- 5.0	ppm
Quan range	+/- 0.5	min
Main isomer peak filter	ON	
ID Quality filter	A, B, C, D	
Adducts (pos. ion)	+H, NH ₄ , Na, +H-H ₂ O	
Adducts (neg. ion)	-H, -CH ₃ , -CH ₂ CO ₂	
Lipid Classes	*Lipids	
Alignment Parameter	Setting	Units
R.T. tolerance	0.10	min
All isomer peak filter	ON	
m-Score threshold	5.0	
ID Quality filter	A, B, C, D	
Configuration	Setting	
Number data points threshold	5	
Intensity baseline, %	0.05	
Intensity ratio threshold	3	
SN ratio threshold	5	

Table 4. Reproducibility of labeled internal standard in biological replicates

Internal Standard (SPLASH)	Calc. m/z	Rt, min	Conc, μg/mL	CV, %
CHE d ₁₈ :1	657.6441	21.8	35.0	6.2
DG 15:0-18:1(d)	587.5506	12.9	1.00	5.7
LPC 18:1(d)	528.3921	3.09	2.50	4.0
LPE 18:1(d)	488.3451	3.21	0.50	5.6
PA 15:0-18:1(d)	667.5169	9.21	0.70	6.6
PC 15:0-18:1(d)	752.6061	9.11	16.0	4.6
PE 15:0-18:1(d)	710.5591	9.55	0.50	4.8
PG 15:0-18:1(d)	741.5537	8.25	3.00	5.7
PI 15:0-18:1(d)	829.5698	7.90	1.00	5.4
PS 15:0-18:1(d)	754.5490	8.07	0.50	4.9
SM d18:1-18:1(d)	737.6397	8.28	3.00	5.3
TG 15:0-18:1(d)-15:0	811.7646	20.6	5.50	4.9

DISCUSSION

WCR larvae contain a high content of triglycerides in addition to phospholipids, sphingolipids, and sterol lipids. In this work, we focused on PC and TG lipids for further MSⁿ characterization. Figure 5 summarizes the information provided by MS³ analysis for 48:1 TG and provides unequivocal evidence for two isomeric TG species that co-eluted under the conditions of the analysis. Characterization of TG mixtures is often not possible using MS² alone since multiple possible combinations fatty acids may exist.

With HCD MS² data, neutral loss of fatty acid and ammonia is the signature fragmentation observed for several classes of neutral lipids including DG, TG, and sterol esters. The Orbitrap ID-X Tribrid mass spectrometer provides an intelligent workflow for monitoring of class-specific product ion or neutral losses and automatically conducting a predefined experiment on the same precursor ion during the instrument cycle. This approach efficiently provides additional information without wasting time scheduling targeted experiments within predefined retention times.

In this study we observed that phenotypical differences related to developmental stage and diet occurred in a wide range of lipids species and concentrations. Figure 6 shows the changes in PI 18:1_18:3 isomer peak areas versus the developmental (PI 18:1_18:3) and diet (PI 18:2/18:2) conditions. A difference in the fatty acyl composition reveals that two distinct phenotypes are observed.

Table 4 summarizes the variation observed for the SPLASH standards spiked into the biological replicates. Very good reproducibility was observed with CVs less than 7% in the extracted ions from the high-resolution MS scan obtained while also performing further lipid characterization using LC-MS dd-MSⁿ.

Table 5 summarizes the number of lipids per sub-class after filtering. The total number of high-quality lipid annotations submitted for further statistical analysis was 972; in 12 positive and 12 negative ion runs a total of 1131 CID MS²/MS³ results were obtained in addition to HCD MS² spectra for 1522 different lipid adduct ions.

CONCLUSIONS

- This LC MSⁿ lipidomics workflow can be applied to any complex biological sample including plasma, plants, tissues, cells, and whole organisms such as insects.
- The Orbitrap ID-X Tribrid mass spectrometer provides a highly sophisticated and customizable workflow for lipidomics and improves the characterization of lipid structures.
- LipidSearch software combines positive and negative ion HCD MS² and CID MS²/MS³ spectra for database searches providing higher confidence in lipid annotation.
- Almost one thousand lipids were annotated from and quantified with good reproducibility.
- Functional lipids such as PI isomers and sterol esters are readily annotated and quantified at low levels. Different phenotypes are observed between isomeric fatty acyl combinations and congeners with varying numbers of double bonds.
- Methods with insufficient structural characterization miss these meaningful biological differences.

REFERENCES

- Increased Depth and Confidence of Lipidome Analysis from Insect Tissues using Chromatography Based Methods with High-resolution Orbitrap MSⁿ. D Gachotte, Y Adelfinskaya, J Gilbert, R Kiyonami, D Peake, Y Yokoi, ThP 544, Proceedings of the 66th ASMS Conference on Mass Spectrometry and Allied Topics, San Diego, California, June 3–7, 2018.
- New Developments in Mass Spectrometry No 4, Tandem Mass Spectrometry of Lipids: Molecular Analysis of Complex Lipids by Robert C Murphy, The Royal Society of Chemistry, Cambridge, UK, 2015.
- Shorthand notation for lipid structures derived from mass spectrometry. G Liebisch, JA Vizcaino, H Köfeler, M Trötzmüller, WJ Griffiths, G Schmitz, F Spener and MJO Wakelam, *J Lipid Res.* 2013, 54, 1523–1530.
- Combined LC/MS platform for analysis of all major stratum corneum lipids, and the profiling of skin substitutes. J. van Smeden, W A Boten, T Hankemeier, R Rissmann, J A Bouwstra, R J Vreeken, *Biochim et Biophys Acta* 2014, 1841, 70–79.

ACKNOWLEDGMENTS

The authors would like to thank Elena Sokol for her expert help in making the LipidSearch database definition changes for sphingolipid nomenclature, as well as her feedback regarding usability, and improving the quality of lipid annotations.

TRADEMARKS/LICENSES

© 2018 Thermo Fisher Scientific Inc. All rights reserved. SPLASH and LipidoMIX are registered trademarks of Avanti Polar Lipids, Inc. LipidSearch is a trademark of Mitsui Knowledge Industry Co., LTD.. All other 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.