

A New Lipid Software Workflow for Processing Orbitrap-based Global Lipidomics Data in Translational and Systems Biology Research

David A Peake¹, Yasuto Yokoi², Junhua Wang¹, Yingying Huang¹, Madalina Oppermann³

Thermo Fisher Scientific, ¹San Jose, USA and ³Stockholm, Sweden; ²Mitsui Knowledge Industry, Tokyo, Japan

Overview

Purpose: We present a new workflow for high-resolution Thermo Scientific™ Orbitrap™-based mass spectrometers for lipidomics using a model system consisting of a wild-type strain vs. knockout for Co-Q production in yeast¹.

Methods: Lipids in yeast mitochondria were analyzed by high resolution LC-MS and MS/MS. Lipid Search® software, an MS² based search using a comprehensive lipid database, was used to identify the lipid species and determine significant differences.

Results: The yeast lipidomics results obtained from the LC/MS data using Lipid Search are comparable to results obtained using infusion lipidomics. We also compared the lipids identified using metabolomics analysis of the same data set – component finding and molecular weight (MW) search for assignment of metabolites and lipids. Due to the complexity of lipid extracts we found that the comprehensive lipid database MS² search method is superior to the accurate mass based MW search for lipidomics.

Introduction

Purpose: We present a new workflow for high-resolution Thermo Scientific™ Orbitrap™-based mass spectrometers for lipidomics using a model system consisting of a wild-type strain vs. knockout for Co-Q production in yeast¹.

Methods: Lipids in yeast mitochondria were analyzed by high resolution LC-MS and MS/MS. Lipid Search® software, an MS² based search using a comprehensive lipid database, was used to identify the lipid species and determine significant differences.

Results: The yeast lipidomics results obtained from the LC/MS data using Lipid Search are comparable to results obtained using infusion lipidomics. We also compared the lipids identified using metabolomics analysis of the same data set – component finding and molecular weight (MW) search for assignment of metabolites and lipids. Due to the complexity of lipid extracts we found that the comprehensive lipid database MS² search method is superior to the accurate mass based MW search for lipidomics.

Methods

Phenotypes of WT (wild-type) and Knockout (KO) Yeast Strains (*S. Cerevisiae*)

WT yeast continue to grow after glucose is exhausted from the media (Diauxic shift point) whereas KO yeast have a defect in Coenzyme Q production and do not grow after the shift. Duplicate biological replicates of WT and KO yeast were collected post shift for metabolomic/lipidomic analyses and analyzed by LC-MS.

Sample Preparation

Yeast were treated with zymolase, homogenized and mitochondria were enriched by differential centrifugation. Mitochondrial protein levels were determined by BCA assay. Mitochondria (~0.25 mg) were extracted 3 times with 400 µL of IPA for 10 min at 4 °C. After centrifugation, supernatants were combined and vacuum dried. Samples were dissolved in 250 µL of 65:35:5 Acetonitrile, Isopropanol, Water with 5 µg/mL 17:0 PG.

Liquid Chromatography–Mass Spectrometry (LC-MS)

Thermo Scientific™ Accela™ 1250 chromatograph and Accela Open autosampler, 10 µL Injection. Column: 2.1 x 100 mm C18, 2.7µm operated at 260 µL/min and 55 °C. The RP HPLC method¹ is described in S. Bird, et al., *Anal. Chem.* **2011**, *83*, 940–949, 6648–6657. A Thermo Scientific™ Q Exactive™ high-resolution Orbitrap mass spectrometer was operated at 70K resolution for electrospray ionization pos. ion LC-MS and 35K for Top5 MS/MS (CE 35).

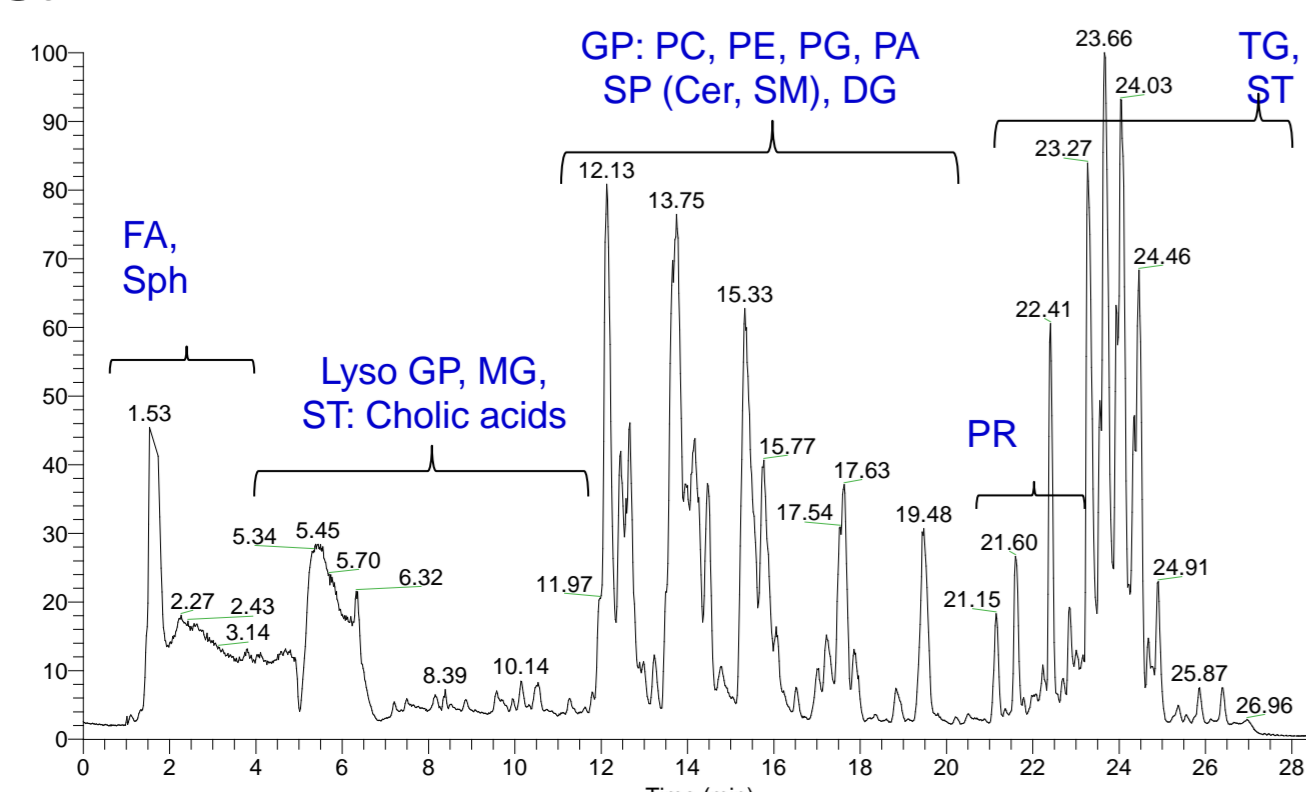
Data Analysis Software

Metabolomics –Thermo Scientific™ SIEVE™ and Lipidomics – Thermo Scientific™ Lipid Search™.

Table 1. Lipid complexity from the LIPID MAPS Structure Database (LMSD)²

Lipid Category	# Class	# Sub-Class	# Lipids
FA	14	36	5,787
GL	6	19	7,568
GP	21	120	6,001
SP	10	31	4,317
ST	6	38	2,678
PR	5	21	1,200
SL	6	7	1,293
PK	15	28	6,741
Total	83	300	37,585

FIGURE 1. LC-MS chromatograms of lipids from WT and KO yeast.

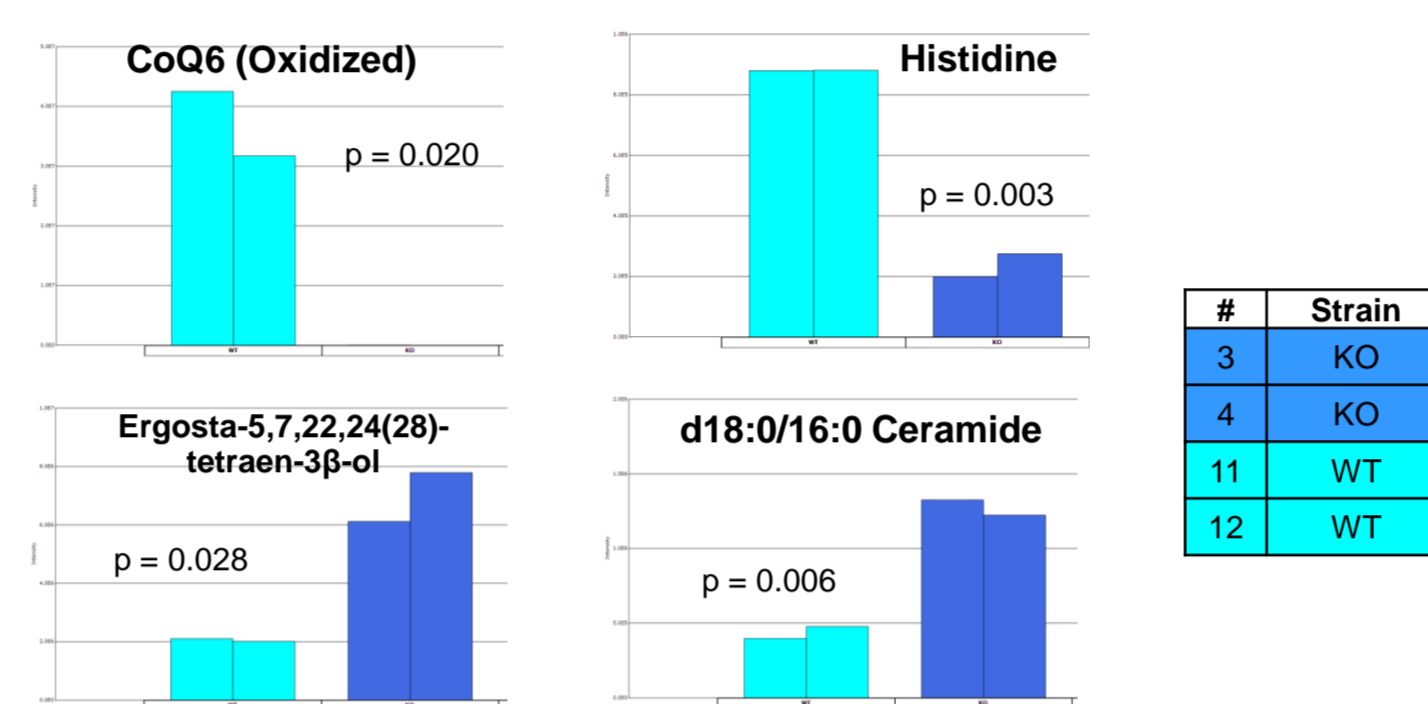


Results

High-Resolution LC-MS Data – Metabolomics Analysis

To characterize the yeast phenotypes we analyzed the sample extracts using an LC-MS method suited for analysis of both metabolites and lipids. The LC-MS chromatogram from WT yeast (Figure 1) shows the regions where lipid classes elute during the LC gradient. Metabolomics analysis using an accurate-mass search tentatively identified 160 metabolites and lipids were present. t-Test statistics (Figure 2) show key metabolite differences.

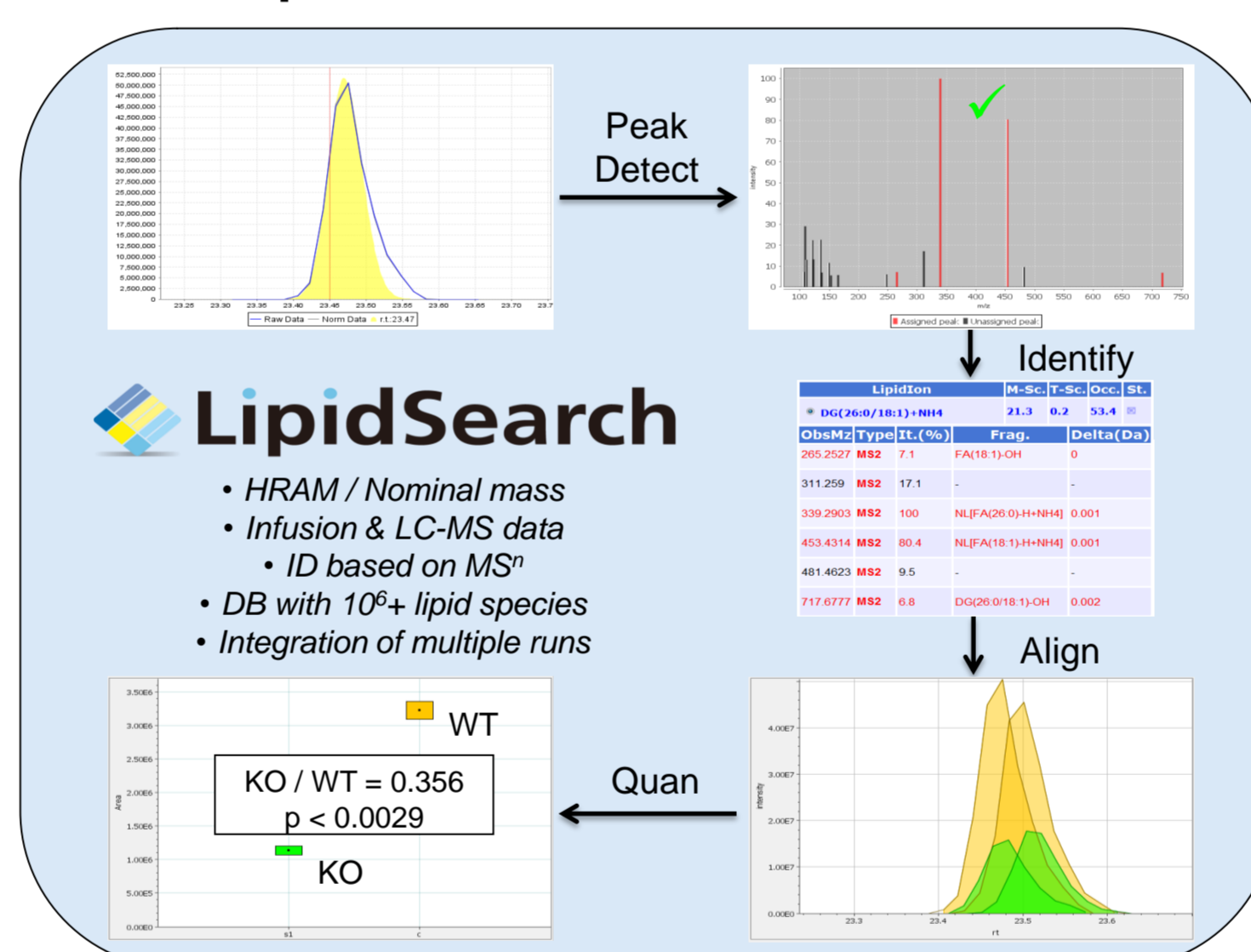
FIGURE 2. Significant metabolite differences observed WT vs. KO yeast.



LC-MS/MS Data Processing Workflow using Lipid Search Software (Figure 3)

- 1) Peak Detection.** Read raw files, MSⁿ and precursor ion accurate masses.
- 2) Identification.** Candidate molecular species are identified by searching a large database > 1,000,000 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 combined 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.
- 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 3. Lipid Search software LC-MS workflow.



Lipid Search Identification and Alignment

LC-MS raw data files containing full scan and data dependent-MS/MS were searched for PL, GL, SP and Co-enzyme lipid classes using a mass tolerance of 5 ppm for precursor ions and 10 ppm for product ions (Figure 4a).

The search results from the 4 samples were aligned using a 0.25 min tolerance window and a combined report was generated (Figure 4b).

FIGURE 4a. Search results for yeast lipids

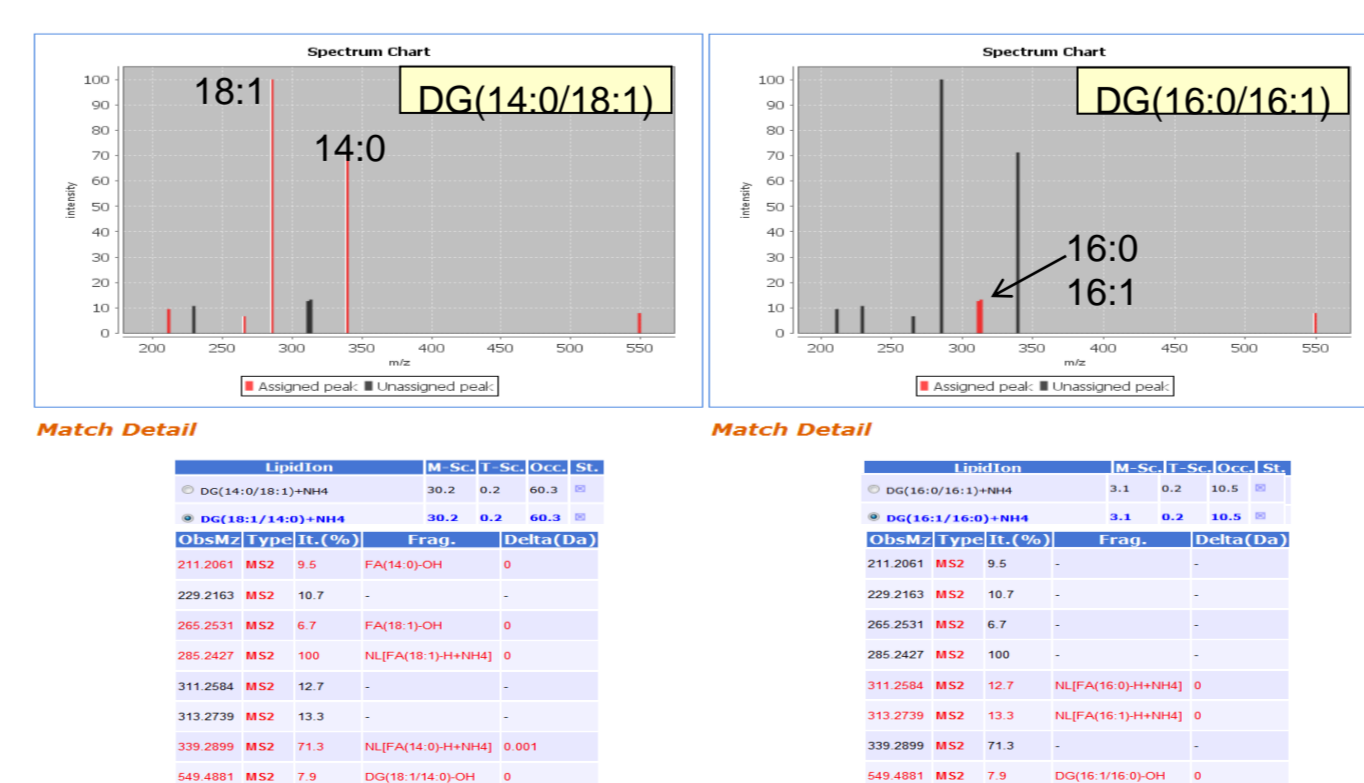
FIGURE 4b. Alignment results for yeast lipids

Search results obtained in < 8 min with 64-bit laptop (MS Windows 7, 2.2 GHz, Intel i7 processor, 8GB RAM)

Identification Report (Figure 5)

For each MS² spectrum, search results are summarized for lipid species matching the predicted fragmentation pattern from the database with a score indicating the fit. If a mixture of lipids is found, the most abundant lipid is displayed. The fragment ions used to identify the lipid are highlighted in red when selected.

FIGURE 5. Search results for m/z 584.5249, Rt = 17.3 min, DG(32:1)



Combined Report – Details (Figures 6 and 7)

Lipid species identified in each LC-MS data file were aligned across the dataset within a retention time tolerance. Quantification is performed on the relative amount of the precursor ion, which in some cases was identified as a mixture of isomers. For each lipid species in the aligned dataset, an interactive report allows review of the data. Relative amounts of each identified lipid were quantified by peak areas and significant differences were determined using t-Tests (Table 2) producing a heat map.

FIGURE 6. Combined report results for PG(17:0/17:0) IS.

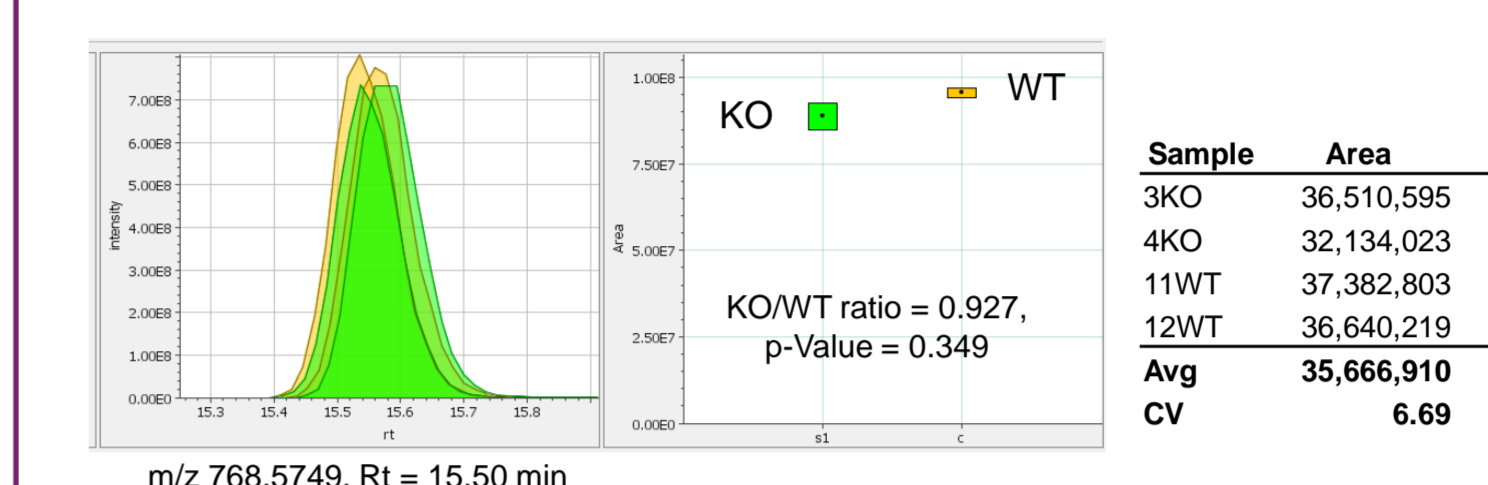
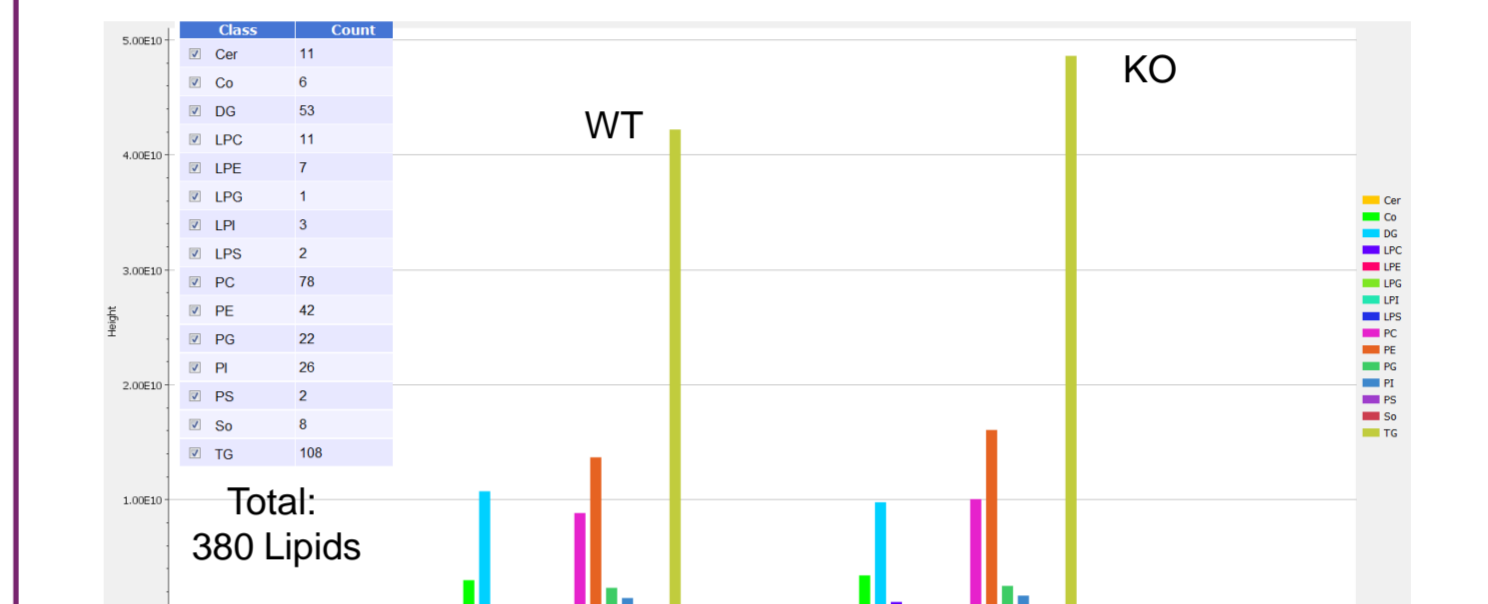


FIGURE 7. Combined report results – total lipid profile.



Yeast Lipidomics Results

The total number of lipids species identified in yeast WT and KO mitochondria (380) is comparable to the number of lipids quantified (250) by infusion lipidomics⁵.

Table 2. Summary of differences between WT vs. KO yeast lipids. Analytes with p-Values < 0.05 for t-Test between WT and KO groups. Fold-change (KO vs. WT) indicated in Red (increase) or Green (decrease).

Class	Compound	RT min	Ratio	p-Value	Class	Compound	RT min	Ratio	p-Value	Class	Compound	RT min	Ratio	p-Value
Cer	Cer(18:0)	15.05	0.006	0.006	CE	CE(18:0)	15.05	0.006	0.006	CE	CE(18:0)	15.05	0.006	0.006
GP	GP(18:0)	15.05	0.006	0.006	GP	GP(18:0)	15.05	0.006	0.006	GP	GP(18:0)	15.05	0.006	0.006
SP	SP(18:0)	15.05	0.006	0.006	SP	SP(18:0)	15.05	0.006	0.006	SP	SP(18:0)	15.05	0.006	0.006
ST	ST(18:0)	15.05	0.006	0.006	ST	ST(18:0)	15.05	0.006	0.006	ST	ST(18:0)	15.05	0.006	0.006
PR	PR(18:0)	15.05	0.006	0.006	PR	PR(18:0)	15.05	0.006	0.006	PR	PR(18:0)	15.05	0.006	0.006
SL	SL(18:0)	15.05	0.006	0.006	SL	SL(18:0)	15.05	0.006	0.006	SL	SL(18:0)	15.05	0.006	0.006
PK	PK(18:0)	15.05	0.006	0.006	PK	PK(18:0)	15.05	0.006	0.006	PK	PK(18:0)	15.05	0.006	0.006

Conclusion

- Lipid Search provides an automated workflow for high quality Orbitrap LC-MS/MS lipidomics data and enables reliable and comprehensive lipid identification.
- Lipid Search identified 380 lipids in MS² spectra from single Orbitrap scans and 112 significant changes were found in the WT and KO yeast phenotypes.
- MS² searching using Lipid Search is a more efficient approach than component finding and MW search for lipid identification.
- Lipid Search reliably identifies product ion mixtures from two or more lipids.
- Data analysis time was dramatically reduced from hours to a few minutes.

References

- Quantitative yeast lipidomics via LC-MS profiling using the Q Exactive Orbitrap mass spectrometer, Peake et al., LIPID MAPS Annual Meeting 2012, May 7-8, 2012, La Jolla, CA.
- LIPID MAPS comprehensive classification system for lipids. E Fahy et al., *J. Lipid Res.* **2009**, *50*, S9-S14.
- Precise and global identification of phospholipid molecular species by an Orbitrap mass spectrometer and automated search engine Lipid Search, R Taguchi, et al., *J. Chrom. A*, **2010**, *1217*, 4229–4239.
- Development of a lipid profiling system using reverse-phase liquid chromatography coupled to high-resolution mass spectrometry with rapid polarity switching and an automated lipid identification software, T Yamada, et al., *J. Chrom. A*, **2013**, *1292*, 211-218.
- Global analysis of the yeast lipidome by quantitative shotgun mass spectrometry, C. Eising, et al., *Proc Natl Acad Sci USA*, **2009**, *102*, 17981–17986.

Acknowledgements

We would like to thank Professor David Pagliarini from the University of Wisconsin, Biochemistry Department for supplying the yeast mitochondrial lipid extracts.

Lipid Search is a registered trademark of MKI, Windows is a trademark of Microsoft and i7 a trademark of Intel. 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. For research use only. Not for use in diagnostic procedures.

