

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

David A Peake,¹ Yasuto Yokoi,² Junhua Wang,¹ and Yingying Huang¹

¹Thermo Fisher Scientific, San Jose, USA; 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

Application of lipidomics to disease phenotype analysis is a growing area in medical research. Identification of unique biomarkers to distinguish healthy humans compared to individuals with disease can have an impact on the early detection of diseases and personalized medicine.

The complexity of the lipidome (Table 1) includes 8 major categories of lipids, over 80 major classes, 300 sub-classes and thousands of lipid species² many with overlapping isomeric or isobaric molecular ions. Because of this complexity, MW searches alone are not sufficient to identify lipids in a complex biological extract.

Identification of lipids requires sophisticated software with an extensive database. The combination of ultra-high resolution MS and MSⁿ analysis should provide unambiguous and precise identification of lipids in biological samples. A robust algorithm for database searching of high-resolution data was developed by Professor Ryo Taguchi and co-workers³ and was commercialized by MKI (Tokyo, JP) as described recently⁴.

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 was 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 (ESI) pos. ion LC-MS and 35K for Top5 MS/MS (CE 35).

Data Analysis Software

Metabolomics –Thermo Scientific™ SIEVE™ software and Lipidomics – Lipid Search software (MKI).

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

	Lipid Category	# Class	# Sub-Class	# Lipids
FA	Fatty acyls	14	36	5,787
GL	Glycerolipids	6	19	7,568
GP	Glycerophospholipids	21	120	8,001
SP	Sphingolipids	10	31	4,317
ST	Sterol lipids	6	38	2,678
PR	Prenol lipids	5	21	1,200
SL	Saccharolipids	6	7	1,293
PK	Polyketides	15	28	6,741
	Total	83	300	37,585

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. Principal component analysis (Figure 2) and t-Test statistics (Figure 3) show key metabolite differences.

FIGURE 1. LC-MS Chromatograms of Lipids from WT and KO Yeast

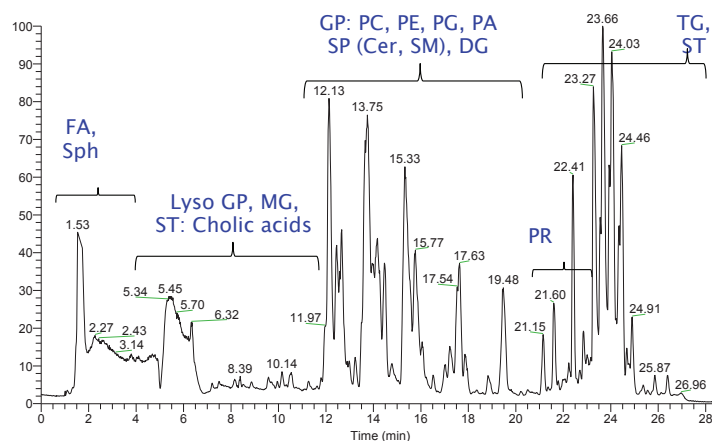


FIGURE 2. Principal Components Analysis of WT and KO Yeast Metabolites

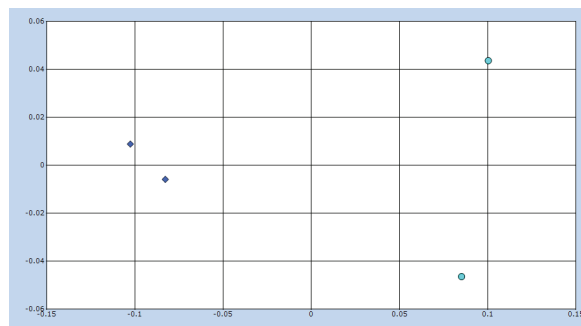
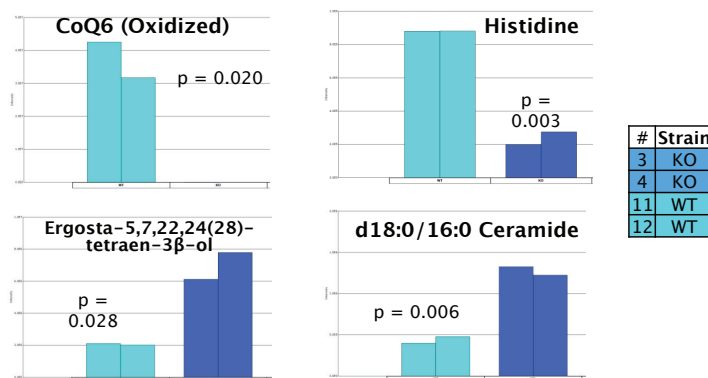


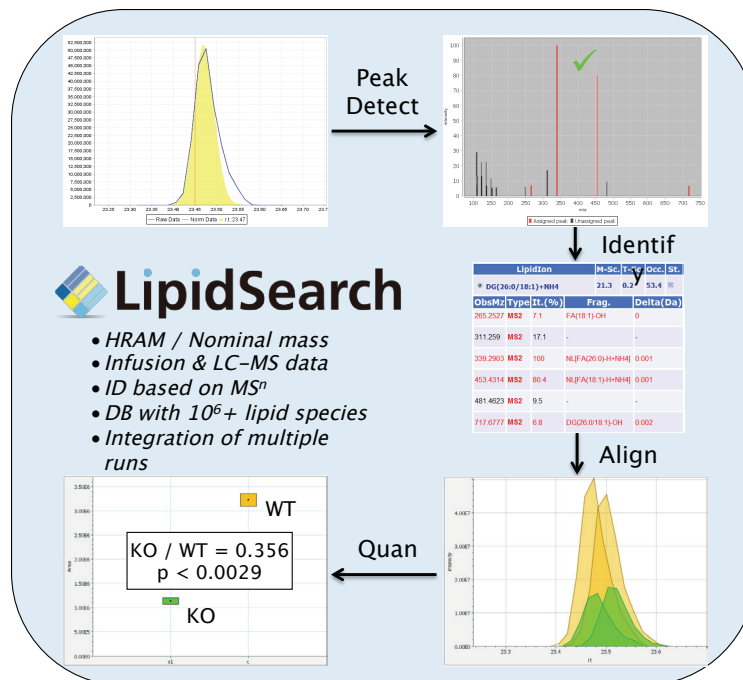
FIGURE 3. Significant Metabolite Differences Observed WT vs. KO Yeast



LC-MS/MS Data Processing Workflow using Lipid Search Software (Figure 4)

- 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 4. Lipid Search Software LC-MS Workflow.



Submitting Data for 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 5a).

The search results from the 4 samples were aligned using a 0.25 min tolerance window and a combined report was generated (Figure 5b).

FIGURE 5a. Search Results for Yeast Lipids

Name	RawData	Type	ExpType	Process	Result	Regist	Update
2013-05-25-21-54-24	4KOPost_1.raw	Product	LC	P I Q	239 %257	2013/05/25 21:54:24	2013/05/25 22:01:01
2013-05-25-21-54-24	3KOPost_1.raw	Product	LC	P I Q	232 %261	2013/05/25 21:54:24	2013/05/25 21:59:38
2013-05-25-21-54-24	12WTPost_1.raw	Product	LC	P I Q	250 %278	2013/05/25 21:54:24	2013/05/25 21:58:12
2013-05-25-21-54-24	11WTPost_1.raw	Product	LC	P I Q	245 %269	2013/05/25 21:54:24	2013/05/25 21:56:58

FIGURE 5b. Alignment Results for Yeast Lipids

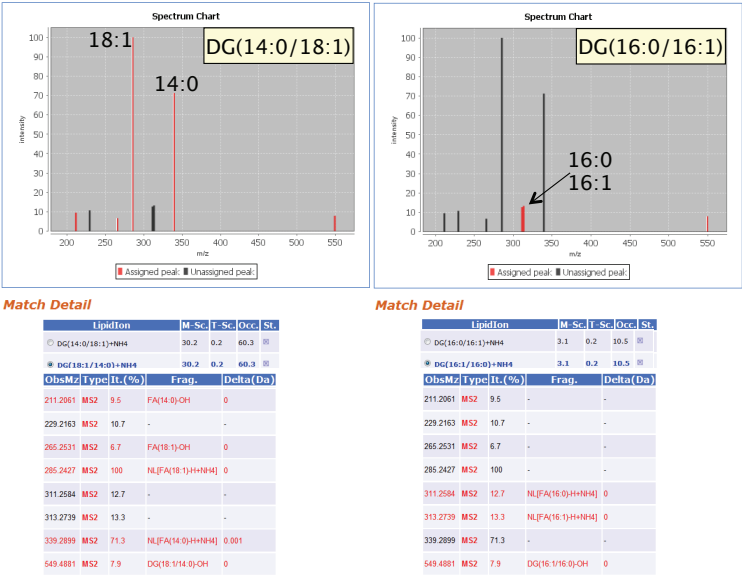
Name	Type	ExpType	Process	Result	Regist	Update
Yeast Lipids	Product	LC	M	380 %700	2013/05/25 22:06:18	2013/05/25 22:07:32

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

Identification Report (Figure 6)

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 each of the species are selected.

FIGURE 6. Search Results for *m/z* 584.5249, *Rt* = 17.3 min, DG(32:1)



Combined Report – Details (Figures 7 and 8)

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 7. Combined Report Results for PG(17:0/17:0) Internal Standard

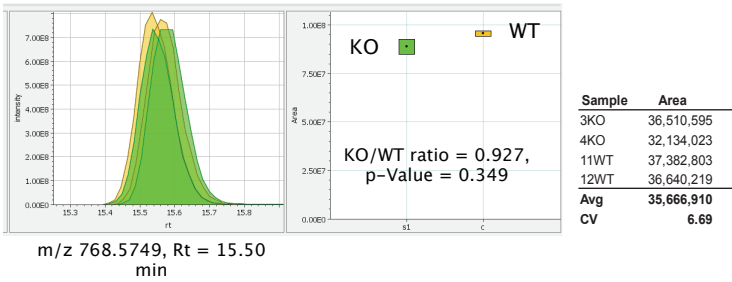
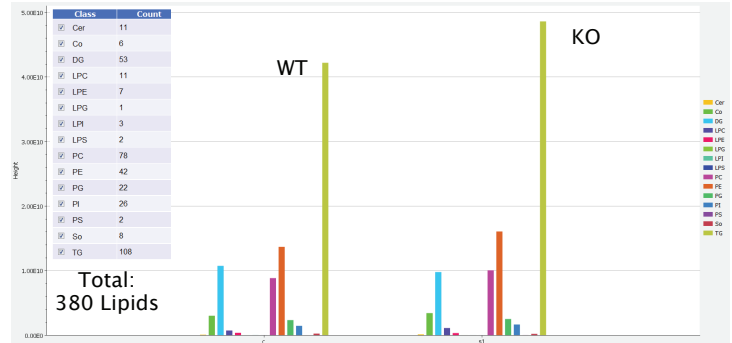


Figure 8. Combined Report Results – Total Lipid Profile



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 by **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(d18:0/16:0)	16.73	2.83	0.006	CoQ	Co(Q6)	15.30	0.09	0.021	DG	DG(16:0/12:0)	15.45	1.33	0.027
	Cer(d18:0/16:1)	15.06	0.52	0.008		Co(Q7)	18.37	0.15	0.017		DG(16:1/15:0)	16.54	0.55	0.027
	Cer(d18:0/18:0)	18.77	2.00	0.037		Co(Q8)	21.15	1.62	0.033		DG(16:1/15:1)	14.84	0.26	0.009
	Cer(d18:0/28:6)	22.74	104.9	0.011		Co(Q9)	22.40	1.50	0.028		DG(16:1/18:3)	14.81	0.53	0.048
	Cer(d18:1/28:6)	22.49	5.93	0.002	PE	PE(10:0/16:0)	10.51	1.83	0.034		DG(16:1/24:0)	22.65	0.35	0.012
	Cer(d18:2/18:1)	16.72	2.07	0.049		PE(10:0/17:1)	10.01	1.65	0.040		DG(18:0/18:1)	21.15	0.45	0.039
So	So(d18:0)	3.03	0.28	0.026		PE(10:0/18:0)	12.55	0.11	0.019		DG(18:1/18:1)	19.54	0.30	0.009
	So(d20:0)	4.79	0.08	0.031		PE(12:0/14:0)	10.51	1.83	0.034		DG(18:1/18:3)	16.60	0.36	0.018
	So(d20:1)	4.97	0.20	0.003		PE(16:0/12:0)	12.25	1.48	0.022		DG(26:0/14:0)	23.10	0.54	0.007
	PC(10:0/16:0)	10.02	6.08	0.000		PE(16:0/15:1)	13.49	1.29	0.018		DG(26:0/16:1)	23.12	0.46	0.023
PC	PC(12:0/18:2)	12.62	2.29	0.004		PE(16:0/16:1)	14.15	1.14	0.028		DG(26:0/18:1)	23.50	0.33	0.005
	PC(12:0/18:2)	12.88	1.50	0.009		PE(16:1/12:0)	10.62	1.83	0.003		DG(26:1/16:1)	22.60	0.06	0.005
	PC(15:0/18:2)	14.36	2.04	0.045		PE(16:1/12:0)	10.96	1.39	0.045		DG(26:1/18:1)	23.05	0.16	0.003
	PC(15:1/12:0)	9.58	2.75	0.023		PE(16:1/15:0)	13.49	1.29	0.018		DG(28:0/18:1)	23.86	0.13	0.000
	PC(16:0/12:0)	11.79	2.95	0.007		PE(16:1/16:1)	12.89	1.57	0.005	TG	TG(10:0/12:0/16:0)	22.25	2.05	0.041
	PC(16:0/12:0)	12.29	3.25	0.024		PE(16:1/16:1)	13.18	1.49	0.021		TG(10:0/14:0/16:0)	22.83	2.52	0.036
	PC(16:0/17:1)	15.71	1.51	0.029		PE(16:1/18:1)	14.26	1.10	0.023		TG(10:0/14:0/16:1)	22.28	3.61	0.042
	PC(16:0/22:6)	12.70	0.23	0.015		PE(17:1/12:0)	11.80	1.54	0.029		TG(10:0/16:0/16:0)	23.28	3.87	0.005
	PC(16:0e/15:1)	18.37	0.11	0.021		PE(18:0/18:2)	15.97	0.25	0.000		TG(10:0/16:0/16:1)	22.84	3.91	0.020
	PC(16:1/12:0)	12.40	3.18	0.003		PE(18:1/14:0)	14.15	1.14	0.028		TG(10:0/16:0/17:1)	23.16	2.44	0.041
	PC(16:1/13:0)	11.31	1.72	0.003		PE(18:1/18:1)	15.91	0.35	0.003		TG(10:0/16:1/16:1)	22.30	3.58	0.032
	PC(16:1/14:0)	14.09	2.08	0.035	PG	PG(16:0/17:1)	13.48	0.95	0.003		TG(12:0/12:0/14:0)	22.25	2.05	0.041
	PC(16:1/16:1)	14.24	1.69	0.002		PG(16:0/18:1)	13.97	0.95	0.007		TG(16:0/12:0/16:0)	23.68	2.08	0.002
	PC(16:1/18:2)	12.71	1.19	0.043		PG(16:0/18:2)	13.08	1.24	0.010		TG(16:0/12:0/16:1)	23.26	2.12	0.013
	PC(16:1/18:3)	11.85	2.47	0.002		PG(16:1/18:1)	12.85	1.28	0.047		TG(16:0/12:0/24:0)	25.39	2.08	0.032
	PC(16:1/20:4)	12.24	0.40	0.007		PG(17:1/17:1)	13.08	1.26	0.010		TG(16:0/14:0/15:0)	23.86	1.37	0.027
	PC(16:1/20:5)	11.12	0.38	0.035		PG(17:1/18:1)	13.53	1.08	0.007		TG(16:0/14:0/16:0)	24.08	1.98	0.004
	PC(17:0/16:0e)	20.60	0.14	0.012	PI	PG(17:1/19:1)	14.63	1.17	0.021		TG(16:0/14:0/16:1)	23.67	1.36	0.029
	PC(17:0/18:0e)	18.37	0.08	0.023		PI(10:0/16:0)	8.56	2.74	0.017		TG(16:0/15:0/16:0)	24.26	1.27	0.022
	PC(18:0/17:1)	17.56	1.56	0.007		PI(12:0/14:0)	8.56	2.74	0.017		TG(16:0/16:0/16:1)	24.07	1.61	0.025
	PC(18:0/18:1)	17.23	0.66	0.045		PI(15:0/18:1)	12.80	0.57	0.043		TG(16:0/16:0/17:0)	24.69	1.21	0.037
	PC(18:0/18:2)	15.48	0.48	0.006		PI(16:1/15:0)	11.23	0.64	0.022		TG(16:0/16:1/16:1)	23.67	1.22	0.046
	PC(18:0/24:2)	21.08	1.85	0.026		PI(16:1/17:0)	12.80	0.57	0.043		TG(16:1/12:0/15:0)	23.16	2.44	0.041
	PC(19:0/18:2)	16.71	0.45	0.003	PS	PI(16:1/18:2)	10.94	0.59	0.047		TG(16:1/18:1/22:1)	24.83	0.62	0.046
	PC(20:0/18:2)	17.20	0.39	0.003		PS(16:1/16:1)	10.86	0.34	0.003		TG(17:1/18:1/18:1)	24.31	0.52	0.017
	PC(20:0/24:1)	22.76	0.51	0.011		PS(16:1/17:1)	11.94	2.26	0.010		TG(18:0/16:0/18:0)	25.39	1.09	0.032
	PC(18:0/18:1)	8.48	2.99	0.015							TG(18:1/18:1/18:1)	24.41	0.50	0.007
	PC(18:0/18:1)	8.80	5.38	0.025							TG(18:1/18:1/18:3)	23.80	0.54	0.019

Conclusions

- 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, D Peake, et al., presented at the 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. doi: 10.1194/jlr.R800095-JLR200.
- 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. doi: 10.1016/j.chroma.2010.04.034.
- 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. doi: 10.1016/j.chroma.2013.01.078
- Global analysis of the yeast lipidome by quantitative shotgun mass spectrometry, C. Ejsing, et al., *Proc Natl Acad Sci USA*, **2009**, *102*, 17981-17986. doi: 10.1073/pnas.0811700106

Acknowledgements

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

www.thermoscientific.com

©2013 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, Inc. and its subsidiaries. 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-Other +27 11 570 1840
Australia +61 3 9757 4300
Austria +43 1 333 50 34 0
Belgium +32 53 73 42 41
Canada +1 800 530 8447
China +86 10 8419 3588
Denmark +45 70 23 62 60

Europe-Other +43 1 333 50 34 0
Finland/Norway/Sweden
+46 8 556 468 00
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9434
Italy +39 02 950 591

Japan +81 45 453 9100
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
Russia/CIS +43 1 333 50 34 0
South Africa +27 11 570 1840



Spain +34 914 845 965
Switzerland +41 61 716 77 00
UK +44 1442 233555
USA +1 800 532 4752

Thermo
SCIENTIFIC
Part of Thermo Fisher Scientific