

# 用于血浆脂质组详细分析的综合、优化、高通量方法

## 作者

Kevin Huynh,  
Natalie A. Mellett,  
Thy Duong, Anh Nguyen,  
Thomas G. Meikle,  
Corey Giles, Peter J. Meikle  
贝克心脏与糖尿病研究所  
澳大利亚墨尔本  
Mark Sartain  
安捷伦科技有限公司

## 摘要

通过将液相色谱/串联质谱 (LC/MS/MS) 等先进脂质组学技术用于群体水平上的人脂质组研究，研究人员得以证实疾病状态与脂类代谢之间存在一定关联。这类研究必然会涉及来自人群的大量样品。全面了解脂质组需要分析数百种不同的脂质，因此需要在覆盖深度和样品通量之间取得平衡。本应用简报介绍了一种基于 LC/TQ 的靶向方法，该方法使用 10  $\mu$ L 血浆测定 763 种不同的人脂质。使用单柱设置时，每个样品的总分析时间为 16 分钟，而运行双柱系统时则缩短至 13 分钟。此外，本文还讨论了样品前处理和脂质萃取的注意事项，以及质量控制策略、内标配制、仪器条件和数据处理。

## 前言

脂质是一类复杂的生物分子，具有多种特性和功能。它们包括细胞膜和脂肪组织的主要成分，并且是哺乳动物系统中的主要能量储存分子。它们还在细胞信号转导通路中发挥着不可或缺的作用。因此，深入了解脂质生物学对于从农业到医学的各个科学领域具有重要意义。

脂质组学是指对脂质组的研究，而脂质组是指给定生物系统（例如细胞、组织或血浆）中的全部脂质。它的最终目标是全面了解存在的脂质种类及其代谢，以及相关生物学信息。近年来，脂质组学技术的发展使我们能够了解脂质组中的各种脂质，包括同分异构体和同质量异构体。特别是液相色谱/串联质谱 (LC/MS/MS) 和多反应监测 (MRM) 等技术，可以帮助用户在短时间内测量数百种不同的脂质，并具有高特异性。新仪器具有更高的灵敏度，有助于测定生物材料中这些浓度极低的化合物。色谱分离性能的提升也进一步增强了更详细地了解脂质组的能力。

对于这些先进的脂质组学技术而言，最令人兴奋的应用是在群体水平研究人类脂质组，进而揭示脂质种类、代谢途径和疾病状态之间的关联。这类研究采用人血浆进行，其中的各种脂质主要存在于脂蛋白颗粒中（图 1），样品来自包含数千名受试者的大型群体和临床研究<sup>[1,2]</sup>。对于如此多的样品，必须在覆盖深度、采集的质谱数据的分辨率以及样品通量之间取得平衡。

考虑到这些因素，本应用简报详细介绍了一种方法，使用 10  $\mu\text{L}$  血浆样品或者等量细胞或组织匀浆测定来自 44 种脂质类别的 763 种不同脂质的浓度。先用反相液相色谱进行分离，然后使用串

联四极杆质谱仪 (MS/MS) 进行分析。关于不同种类/类别脂质的离子化和碎裂的详细信息以及所使用的内标详见表 1。使用单色谱柱方法时，每个样品的总分析时间为 16 分钟，而使用双柱系统的运行时间则为 13 分钟。本文中的方法使用与 Agilent MassHunter 软件包集成的安捷伦 LC/MS/MS 仪器，可实现快速数据处理。

虽然所报告的方法主要为大型队列研究实现高通量血浆样品分析而开发，但通过简单调整后也可用于各种生物基质。也可以根据所需的结构分辨率调整 LC 分离和 MS/MS 参数，或根据通量要求修改分离方案。

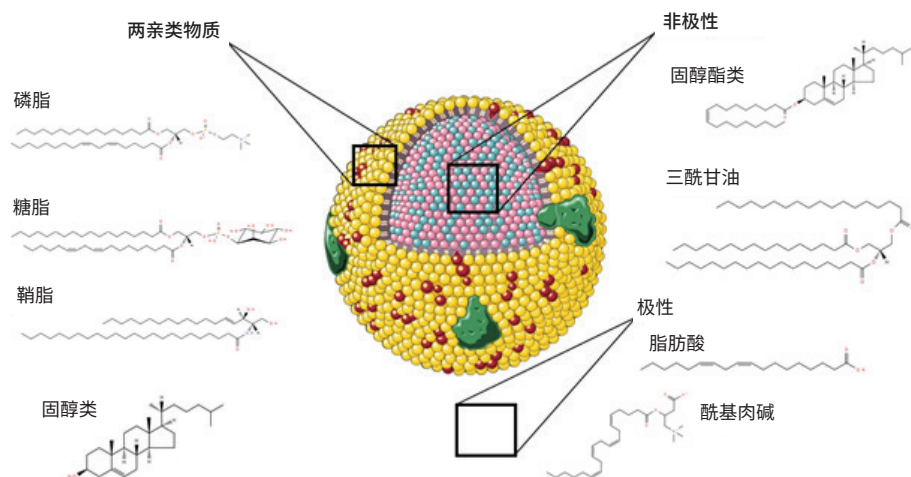


图 1. 人血浆中常见脂质类别的示意图。脂质是一类小型两亲分子，包含多种类别和亚类。在人血浆中，这些脂质主要存在于外周循环的脂蛋白颗粒中

表 1. 不同脂质类别离子化和碎裂的详细信息

脂质类别/亚类	母离子	碎裂 <sup>1</sup>	脂质种类数量	内标	内标 (pmol)/样品
鞘氨醇 (Sph)	[M + H] <sup>+</sup>	NL, 18.0 Da	2	Sph(d17:1)	20
鞘氨醇-1-磷酸 (S1P)	[M + H] <sup>+</sup>	特定鞘氨醇碱	4	S1P(18:1) d7	40
二氢神经酰胺 (dhCer)	[M + H] <sup>+</sup>	特定鞘氨醇碱	6	dhCer(d18:0/8:0)	50
神经酰胺 (Cer(d))	[M + H] <sup>+</sup>	特定鞘氨醇碱	41	Cer(d18:1-d7/18:0)	50
脱氧神经酰胺 (Cer(m))	[M + H] <sup>+</sup>	特定鞘氨醇碱	11	Cer(d18:1-d7/18:0)	50
神经酰胺-1-磷酸酯 (C1P)	[M + H] <sup>+</sup>	特定鞘氨醇碱	1	Cer(d18:1-d7/18:0)	50
单己糖神经酰胺 (HexCer)	[M + H] <sup>+</sup>	特定鞘氨醇碱	13	HexCer(d18:1/15:0) d7	50
双己糖神经酰胺 (Hex2Cer)	[M + H] <sup>+</sup>	特定鞘氨醇碱	9	Hex2Cer(d18:1/15:0) d7	50
三己糖神经酰胺 (Hex3Cer)	[M + H] <sup>+</sup>	特定鞘氨醇碱	5	Hex3Cer(d18:1/17:0)	50
GM3 神经节苷脂 (GM3)	[M + H] <sup>+</sup>	特定鞘氨醇碱	6	Hex3Cer(d18:1/17:0)	50
硫苷脂 (SHexCer)	[M + H] <sup>+</sup>	PI, m/z 264.3	5	SHexCer(d18:1/12:0)	10
鞘磷脂 (SM)	[M + H] <sup>+</sup>	PI, m/z 184.1	45	SM(d18:1/15:0) d9	100
磷脂酸 (PA)	[M + H] <sup>+</sup>	NL, 115.0 Da	4	PA(15:0_18:1) d7	50
磷脂酰胆碱 (PC)	[M + H] <sup>+</sup>	PI, m/z 184.1	70	PC(15:0_18:1) d7	100
烷基磷脂酰胆碱 (PC(O))	[M + H] <sup>+</sup>	PI, m/z 184.1	23	PC(15:0_18:1) d7	100
烯基磷脂酰胆碱 (缩醛磷脂) (PC(P))	[M + H] <sup>+</sup>	PI, m/z 184.1	28	PC(P-18:0/18:1) d9	100
溶血磷脂酰胆碱 (LPC)	[M + H] <sup>+</sup>	PI, m/z 184.1 和 m/z 104.1	61	LPC(18:1) d7	100
溶血烷基磷脂酰胆碱 (溶血血小板活化因子) (LPC(O))	[M + H] <sup>+</sup>	PI, m/z 104.1	10	LPC(18:1) d7	100
溶血烯基磷脂酰胆碱 (缩醛磷脂) (LPC(P))	[M + H] <sup>+</sup>	PI, m/z 104.1	6	LPC(18:1) d7	100
磷脂酰乙醇胺 (PE)	[M + H] <sup>+</sup>	NL, 141.0 Da	32	PE(15:0_18:1) d7	100
烷基磷脂酰乙醇胺 (PE(O))	[M + H] <sup>+</sup>	NL, 141.0 Da	15	PE(15:0_18:1) d7	100
烯基磷脂酰乙醇胺 (缩醛磷脂) (PE(P))	[M + H] <sup>+</sup>	特定酰基链	48	PE(P-18:0/18:1) d9	100
溶血磷脂酰乙醇胺 (LPE)	[M + H] <sup>+</sup>	NL, 141.0 Da	12	LPE(18:1) d7	100
溶血烯基磷脂酰乙醇胺 (缩醛磷脂) (LPE(P))	[M + H] <sup>+</sup>	NL, 171.9 Da	4	LPE(18:1) d7	100
磷脂酰肌醇 (PI)	[M + NH <sub>4</sub> ] <sup>+</sup>	NL, 277.0 Da	25	PI(15:0_18:1) d7	50
磷脂酰肌醇单磷酸 (PIP1)	[M + NH <sub>4</sub> ] <sup>+</sup>	NL, 357.0 Da	1	PI(15:0_18:1) d7	50
溶血磷脂酰肌醇 (LPI)	[M + NH <sub>4</sub> ] <sup>+</sup>	NL, 277.0 Da	7	LPI(13:0)	20
磷脂酰丝氨酸 (PS)	[M + H] <sup>+</sup>	NL, 185.0 Da	4	PS(15:0_18:1) d7	50
磷脂酰甘油 (PG)	[M + NH <sub>4</sub> ] <sup>+</sup>	NL, 189.0 Da	4	PG(15:0_18:1) d7	50
胆固醇酯 (CE)	[M + NH <sub>4</sub> ] <sup>+</sup>	PI, m/z 369.3	27	CE(18:0) d6	1000
游离胆固醇 (COH)	[M + NH <sub>4</sub> ] <sup>+</sup> 源内碎片	PI, m/z 161.2	1	COH(d7)	10000
脱氢胆固醇酯 (DE)	[M + NH <sub>4</sub> ] <sup>+</sup>	PI, m/z 367.4	4	CE(18:0) d6	1000
甲基-胆固醇酯 (甲基-CE)	[M + NH <sub>4</sub> ] <sup>+</sup>	PI, m/z 383.3	4	CE(18:0) d6	1000
甲基-脱氢胆固醇酯 (甲基-DE)	[M + NH <sub>4</sub> ] <sup>+</sup>	PI, m/z 381.4	2	CE(18:0) d6	1000
二甲基-胆固醇酯 (二甲基-CE)	[M + NH <sub>4</sub> ] <sup>+</sup>	PI, m/z 396.6	3	CE(18:0) d6	1000
游离脂肪酸 (FFA)	[M-H] <sup>-</sup>	SIM	16	FFA(18:1) d9	200
酰基肉碱 (AC)	[M + H] <sup>+</sup>	PI, m/z 85.1	28	AC(16:0) d3	10
羟基化酰基肉碱 (AC-OH)	[M + H] <sup>+</sup>	PI, m/z 85.1	8	AC(16:0) d3	10
胆汁酸 (BA)	[M + NH <sub>4</sub> ] <sup>+</sup>	取决于种类	2	胆酸 d4	50
二酰甘油 (DG)	[M + NH <sub>4</sub> ] <sup>+</sup>	NL, 脂肪酸	24	DG(15:0_18:1) d7	200
三酰甘油 (中性丢失, 关联研究) (TG [NL])	[M + NH <sub>4</sub> ] <sup>+</sup>	NL, 脂肪酸	77	TG(15:0_18:1-d7_15:0)	100
烷基二酰甘油 (中性丢失, 关联研究) (TG(O) [NL])	[M + NH <sub>4</sub> ] <sup>+</sup>	NL, 脂肪酸	19	TG(15:0_18:1-d7_15:0)	100
泛醌	[M + NH <sub>4</sub> ] <sup>+</sup>	PI, m/z 197.0	1	Hex3Cer(d18:1/17:0)	50
氧化脂质 (氧化类)	取决于种类 [+O]	取决于种类	6	取决于种类	-

<sup>1</sup> NL = 中性丢失, PI = 母离子

## 实验部分

### 化学品

甲酸铵购自 Sigma-Aldrich (St. Louis, MO, USA)。甲醇 (LC 级)、1-丁醇 (LC 级)、水 (LC/MS 级)、乙腈和异丙醇 (hypergrade, LC/MS 级, LiChrosolv) 均购自 Supelco (Bellefonte, PA, USA)。NIST1950 QC 物质购自美国国家标准与技术研究院 (Los Angeles, CA, USA)。除酰基肉碱 16:0 d3 (Larodan AB, Solna, Sweden)、胆固醇酯 18:0 d6 (CDN

Isotopes, Quebec, Canada)、C18:1 d9 脂肪酸 (Sigma-Aldrich, St. Louis, MI, USA) 和三己糖神经酰胺 17:0 (Matreya LLC, State College, PA, USA) 外, 所有内标均购自 Avanti Polar Lipids (Alabaster, AL, USA)。

### 制备内标混合物

为了测定所分析的每种脂质的相对浓度, 将内标加入每个样品中, 包括稳定同位素标记的内标和非生理性脂质内标。使用市售材料自行制备内标混合物 (表 2), 但预

配制的混合物 (例如 SPLASH LIPIDOMIX 质谱标准品 (Avanti)) 同样适用。如果要使用制备的其他内标, 则需要生成每个新增标准化合物的保留时间, 并将其用于更新仪器和数据处理方法文件。

从供应商获得的各种内标均在  $-80^{\circ}\text{C}$  下储存。配制混标时, 将标准品解冻至室温, 然后复溶于 1:1 氯仿:甲醇中, 最终获得如表 2 第 3 列中所示的规定储备液浓度。然后将所得储备液混合, 得到一份含所有内标的混合物, 各化合物的最终浓度

表 2. 内标信息

名称	分子量	储备液 (μmol/L)	混标中的浓度 (μmol/L)	pmol/样品 (10 μL)	溶剂	供应商	货号	Avanti SPLASH LIPIDOMIX <sup>1,2,3</sup> 中是否存在
AC(16:0) d3	439.09	100	1	10	CHCl <sub>3</sub> :MeOH (1:1)	Larodan	71-1746-5	否
CE(18:0) d6	659.16	2000	100	1000	CHCl <sub>3</sub> :MeOH (1:1)	CDN isotopes	D-5823	否
Cer(d18:1-d7/18:0)	572.99	200	5	50	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	860677P	否
胆酸 d4	412.61	200	5	50	CHCl <sub>3</sub> :MeOH (1:1)	Cayman	20849	否
COH(d7)	393.69	20000	1000	10000	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	700041P	是
DG(15:0_18:1) d7	587.97	800	20	200	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	791647	是
dhCer(d18:0/8:0)	427.71	200	5	50	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	860626P	否
FFA(18:1) d9	291.52	3000	20	200	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	861809	否
Hex2Cer(d18:1/15:0) d7	855.21	400	5	50	CHCl <sub>3</sub> :MeOH:H <sub>2</sub> O (1:1:0.2)	Avanti	330727	否
Hex3Cer(d18:1/17:0)	1038.35	100	5	50	CHCl <sub>3</sub> :MeOH:H <sub>2</sub> O (1:1:0.2)	Matreya LLC	1523	否
HexCer(d18:1/15:0) d7	693.06	400	5	50	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	330729	否
LPC(18:1) d7	528.71	400	10	100	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	791643	是
LPE(18:1) d7	486.63	400	10	100	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	791644	是
LPI(13:0)	547.57	200	2	20	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	850101	否
PA(15:0_18:1) d7	689.93	400	5	50	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	791642	是
PC(15:0_18:1) d7	753.09	400	10	100	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	791637	是
PC(P-18:0/18:1) d9	781.19	400	10	100	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	852475	是
PE(15:0_18:1) d7	711.01	400	10	100	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	791638	是
PE(P-18:0/18:1) d9	739.11	400	10	100	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	852474	是
PG(15:0_18:1) d7	764.01	400	5	50	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	791640	是
PI(15:0_18:1) d7	847.12	400	5	50	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	791641	是
PS(15:0_18:1) d7	777.00	400	5	50	CHCl <sub>3</sub> :MeOH:H <sub>2</sub> O (1:1:0.2)	Avanti	791639	是
S1P(18:1) d7	386.52	400	4	40	CHCl <sub>3</sub> :MeOH:H <sub>2</sub> O (1:1:0.2)	Avanti	860659	否
SHexCer(d18:1/12:0)	741.03	200	1	10	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	860573P	否
SM(d18:1/15:0) d9	698.06	800	10	100	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	860686	否
Sph(d17:1)	285.47	400	2	20	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	860640P	否
TG(15:0_18:1-d7_15:0)	812.35	400	10	100	CHCl <sub>3</sub> :MeOH (1:1)	Avanti	791648	是

<sup>1</sup> Avanti SPLASH LIPIDOMIX 中不包含此处所列的特定 CE 和 SM 标准品, 但包含等效标准品

<sup>2</sup> 缩醛磷脂标准品仅以 SPLASH II Lipidomix 混标的形式提供。胆固醇 (COH)、PA 和 PG 标准品仅以 SPLASH I LIPIDOMIX 的形式提供

<sup>3</sup> SPLASH LIPIDOMIX 中的所有标准品并非都具有第 3 列中所示的浓度

列于表 2 第 4 列。将该内标混合物置于超声仪水浴中超声处理 15 分钟，确保完全混合，然后进行分装以便单次使用，并在 -80 °C 下储存。

### 制备萃取溶剂

为制备用于血浆样品的脂质萃取溶剂，将上述分装的内标混合物解冻至室温（1 小时），超声水浴处理 15 分钟，然后涡旋混合。然后在真空中干燥所需体积的内标混合物（每个待萃取样品 10 µL），并将其复溶于含有 10 mmol/L 甲酸铵的十倍体积的 1:1 丁醇:甲醇中，再次超声处理 15 分钟。例如，如果要萃取 100 个样品，则干燥 1 mL 内标混合物，并将其复溶于含有 10 mmol/L 甲酸铵的 10 mL 1:1 丁醇:甲醇中。

### 样品任务列表、随机分配以及加入质控样品

在萃取或分析样品之前生成了一个样品任务列表，其中指定了试样和对照样品的处理顺序。样品保持这个顺序，首先进行萃取，然后将其引入 LC/MS/MS 进行分析。为了实现这一目的，最初对所有试样进行随机分配，尽可能降低整个数据集的萃取偏差影响。具有统计相关性的样品组或时间点保持配对，而采集顺序等因素则随机分配。

随后，将多个不同的质控样品加入任务列表中。PQCs、NIST1950 QCs、空白样品和试样一起使用相同的程序萃取，而 TQCs 则从之前提取并合并的血浆中提取。这些样品需要比对结果数据，并补偿仪器漂移和误差（批次内和批次间），以及检测数据中更严重的错误。下文总结了每种质量控制类型的组成和作用，以及将它们分散在样品任务列表中所用的程序。

### 血浆质量控制 (PQC)

PQC 样品由采集的健康志愿者 (n = 8-20) 血浆组成，将它们合并后再分装以便单次使用，并在 -80 °C 下储存。为了评估萃取过程中引入的任何误差，每隔 15 个样品加入 1 个或多个 PQC 样品（每个批次至少包含 10 个 PQC 样品）。通过这些样品的 %CV 可以监测移液和处理过程中引入的误差，以及在大量样品需要进行多批次萃取时进行批次比对。

### 技术质量控制 (TQC)

通过萃取用于 PQC 的相同血浆的等分试样，然后合并提取物以消除萃取差异，制得 TQC 样品。在消除萃取差异后，这些 TQC 样品可实现对仪器差异和性能（分离度、保留时间漂移等）的特定监测。每隔 15 个生物样品加入 1 个或多个 TQC 样品，每个批次至少包含 10 个 TQC 样品。

### NIST1950 质量控制 (NIST1950 QC)

NIST1950 是一种市售标准化血浆来源。添加 NIST1950 质控样品不仅有助于准确比对实验室内的不同样品组，还有助于比对同样使用 NIST1950 QC 的其他实验室的不同样品组。为避免多次冻融循环，将 NIST1950 QC 样品解冻并等分 (10 µL) 装入 Eppendorf 萃取管中。每隔 30 个样品加入 1 个或多个 NIST1950 QC 样品，每个批次至少包含 8 个 NIST1950 QC 样品。

### 空白样品

空白样品由 10 µL milliQ 水组成，加入空白样品是为了监测背景信号，空白样品与生物样品和 PQC 同时进行萃取。每隔 30 个样品加入 1 个或多个空白样品，每个批次至少包含 5 个空白样品。

### 人血浆脂质提取

为防止高浓度的盐和蛋白质对仪器造成污染，将所有血浆样品提取到溶剂中，使目标化合物转移到溶液中，而盐和蛋白质则可以沉淀并丢弃<sup>[9]</sup>。

使用聚丙烯外置活塞式移液器在 1.5 mL 微量离心管中进行脂质提取。将 10 µL 血浆与 100 µL 萃取溶剂混合，萃取溶剂由含有 10 mmol/L 甲酸铵的丁醇:甲醇 (1:1) 和复溶的内标混合物组成（表 2）。然后将每个样品涡旋混合 5 秒，并超声水浴处理 1 小时，温度保持在 21-25 °C。然后

对样品进行离心 (13000 xg, 10 分钟, 20 °C), 并将上清液转移至带有 200 µL 玻璃内插管 (5183-2085) 的 1.5 mL 玻璃样品瓶 (5190-9062) 中。应使用 PTFE/硅橡胶盖 (5185-5820) 将样品密封。样品在 -80 °C 下储存, 在分析前才能将其取出, 并立即解冻至室温 (1 小时), 超声水浴处理 15 分钟, 短暂涡旋混合, 然后加载到仪器的自动进样器中。

如果需要高容量分析, 可以按照应用简报 5991-5724<sup>[4]</sup> 所述, 在 Bravo 平台上使用 96 孔板进行样品萃取操作。

### 仪器和 LC/MS/MS 方法

使用 Agilent 1290 Infinity/Infinity II 液相色谱系统分离提取的血浆样品, 该系统包括:

- Agilent 1290 Infinity 自动进样器 (G4226A) **注:** 也可以使用对保留时间影响更小的 1290 Infinity II Multisampler (G7167B)
- Agilent 1290 Infinity II 高速 (二元) 泵 (G7120A, 双柱设置使用两台泵)
- Agilent 1290 Infinity II 大容量柱温箱 (G7116B), 配备 Agilent InfinityLab Quick Change 2 位/10 通快速切换阀头 (用于双柱设置, 部件号 5067-4240)

单柱和双柱设置的管路要求如图 2 和表 5 所示。

该液相色谱系统与配备安捷伦喷射流技术离子源的 Agilent 6495C 三重四极杆液质联用系统联用。

脂质提取物 (1 µL, 相当于 0.1 µL 血浆) 进样, 并采用 16 分钟梯度进行分离, 分离条件详见表 3, 其中前 0.8 分钟导入废液瓶。也可以使用双柱设置, 通过 13 分钟梯度程序 (详见表 4) 来缩短分析时

表 3. 单柱设置的色谱条件

参数	Agilent 1290 Infinity/Infinity II LC																						
分析柱	Agilent ZORBAX Eclipse Plus C18, 100 × 2.1 mm, 1.8 µm (部件号 959758-902)																						
在线过滤器	Agilent 1290 Infinity II 在线过滤器, 0.3 µm (部件号 5067-6189)																						
柱温	45 °C																						
进样量	1 µL (相当于 0.1 µL 血浆)																						
自动进样器温度	20 °C																						
进样针清洗	清洗瓶, 3 s, 1:1 丁醇:甲醇 * 也可以使用清洗口, 10 s, 1:1 流动相 A/B																						
流动相	A) 10 mmol/L 甲酸铵, 5 µmol/L 安捷伦去活剂添加剂 (部件号 5191-3940) 的 5:3:2 水:乙腈:异丙醇溶液 B) 10 mmol/L 甲酸铵的 1:9:90 水:乙腈:异丙醇溶液 * 请勿向流动相 B 中添加去活剂添加剂																						
流速	0.4 mL/min																						
梯度程序	<table border="1"> <thead> <tr> <th>时间</th> <th>%B</th> </tr> </thead> <tbody> <tr><td>0</td><td>15</td></tr> <tr><td>2.5</td><td>50</td></tr> <tr><td>2.6</td><td>57</td></tr> <tr><td>9</td><td>70</td></tr> <tr><td>9.1</td><td>93</td></tr> <tr><td>11</td><td>96</td></tr> <tr><td>11.1</td><td>100</td></tr> <tr><td>12</td><td>100</td></tr> <tr><td>12.2</td><td>15</td></tr> <tr><td>16</td><td>15</td></tr> </tbody> </table>	时间	%B	0	15	2.5	50	2.6	57	9	70	9.1	93	11	96	11.1	100	12	100	12.2	15	16	15
时间	%B																						
0	15																						
2.5	50																						
2.6	57																						
9	70																						
9.1	93																						
11	96																						
11.1	100																						
12	100																						
12.2	15																						
16	15																						
停止时间	16 min																						
后运行时间	无																						

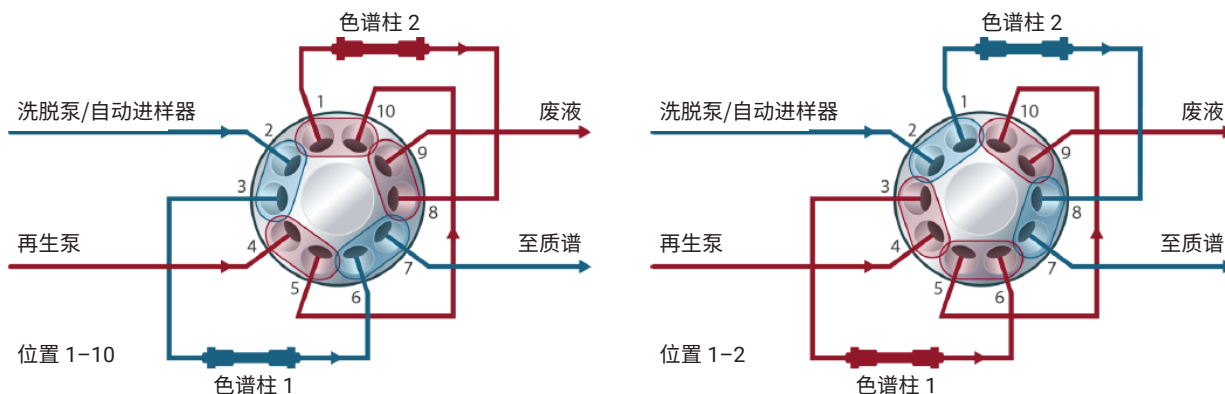


图 2. 双柱设置。通过切换阀管路实现交替色谱柱再生的示例

间。采用动态多反应监测 (dMRM) 进行 LC/MS/MS 分析。MS 仪器参数和离子源条件如表 6 所示。有关该方法的开发及更多详细信息，请参阅之前发表的文章<sup>[1]</sup>。

表 4. 双柱设置的色谱条件

参数	Agilent 1290 Infinity/Infinity II LC		
分析柱	两根 Agilent ZORBAX Eclipse Plus C18, 100 × 2.1 mm, 1.8 μm (部件号 959758-902)		
在线过滤器	两个 Agilent 1290 Infinity II 在线过滤器, 0.3 μm (部件号 5067-6189)		
柱温	45 °C		
进样量	1 μL (相当于 0.1 μL 血浆)		
自动进样器温度	20 °C		
进样针清洗	清洗瓶, 3 s, 1:1 丁醇:甲醇 * 也可以使用清洗口, 10 s, 1:1 流动相 A/B		
流动相	A) 10 mmol/L 甲酸铵, 5 μmol/L 安捷伦去活剂添加剂 (部件号 5191-3940) 的 5:3:2 水:乙腈:异丙醇溶液 B) 10 mmol/L 甲酸铵的 1:9:90 水:乙腈:异丙醇溶液 * 请勿向流动相 B 中添加去活剂添加剂		
流速	0.4 mL/min		
梯度程序	泵 A (G7120A)		
	时间	%B	
	0	15	
	2.5	50	
	2.6	57	
	9	70	
	9.1	93	
	11	96	
	11.1	100	
	12	100	
梯度程序	泵 B (G7120A)		
	时间	%B	
	0	15	
	1	15	
	2	100	
	6	100	
	7	15	
	12.9	15	
	停止时间	13 min	
	后运行时间	无	
MCT 阀位置	使用当前位置, 运行后增加阀位		

表 5. 单柱设置的液相色谱管线详细信息

位置	部件描述	部件号 <sup>1</sup>
二元泵混合器	V35 Jet Weaver 混合器	G7120-68135
从泵到自动进样器的管线	不锈钢毛细管, 内径 0.17 mm, 600 mm	G1312-67305
针头	针头组件, 0.12 mm	G4226-87201
针座	针座, 内径 0.12 mm	G4226-87012
流量计毛细管	不锈钢毛细管, 0.17 × 160 mm, S/SL, ps/ps	G4226-60301
样品定量环组件	20 μL 定量环组件	G4226-60310
自动进样器到 MCT	不锈钢毛细管, 0.12 × 500 mm, 带长套管	5500-1157
热交换器	标准流速热交换器, 1.6 μL	G7116-60015
连接到在线过滤器	Quick Connect 快速连接组件, 不锈钢, 0.12 × 105 mm	5067-5957
热交换器进/出口接头	Quick Turn 液相色谱接头	5067-5966
PEEK 管线到 MS	PEEK 管线, 内径 0.13 mm, 700 mm	0890-1915

<sup>1</sup> 部件号特定用于此处所述的液相色谱配置。如果使用 Agilent 1290 Infinity II Multisampler, 某些部件号会有所不同

## 动态多反应监测设置

该方法的一个关键是在质谱仪中使用动态多反应监测 (dMRM)。如果仪器要测量广泛的脂质组中的大量不同脂质，必须定义每种化合物洗脱的特定保留时间窗口，使仪器仅在整个样品运行时间内的特定点扫描给定的脂质种类。该保留时间窗口因不同的 MRM 而异，并且主要取决于峰宽的差异。这种方法可改善仪器的占空比，并延长每种化合物的平均驻留时间 (图 3)。

使用 Agilent MassHunter 采集软件准备 dMRM 方法，其中每种目标化合物的 MRM 扫描与化合物的已知保留时间大致匹配。保留时间见表 A1 (参见附录)。使用每个脂质类别的代表性标准品优化碰撞能量和 MRM 离子对，尽可能提高响应。使用 Agilent Source Optimizer 软件优化 AJS 离子源参数，平衡所有脂质类别的灵敏度。选择 650 ms 作为循环时间，使整个峰中保持足够数量的数据点，从而实现可靠的定量。本研究利用 665 个

表 6. MS 仪器参数

参数	Agilent 6495C 三重四极杆液质联用系统
离子源	安捷伦喷射流
极性	正负切换
气体温度	150 °C
干燥气 (氮气)	17 L/min
雾化器气体	20 psi
鞘气	200 °C
鞘气流速	10 L/min
毛细管电压	3500 V (+), -3000 V (-)
喷嘴电压	1000 V (+), -1500 V (-)
iFunnel 高/低压力 RF	200/110 V (+), 150/60 V (-)
扫描类型	动态 MRM (dMRM)
Q1/Q2 分辨率	单位 (0.7 amu)
Delta EMV	100 V (+), 0 V (-)
碰撞池加速电压	4-5 V
循环时间	650 ms
时间过滤器宽度	0.07 min
MRM 总数	665 (正离子: 646, 负离子: 19)

离子对来监测 763 种人脂质和异构体。在 665 个离子对中，最“繁忙”的部分得到了 115 个离子对，平均驻留时间不

到 5 ms。即使驻留时间很短，6495C LC/TQ 仍然能够在这些化合物中保持出色的精密度 (参见结果与讨论部分)。

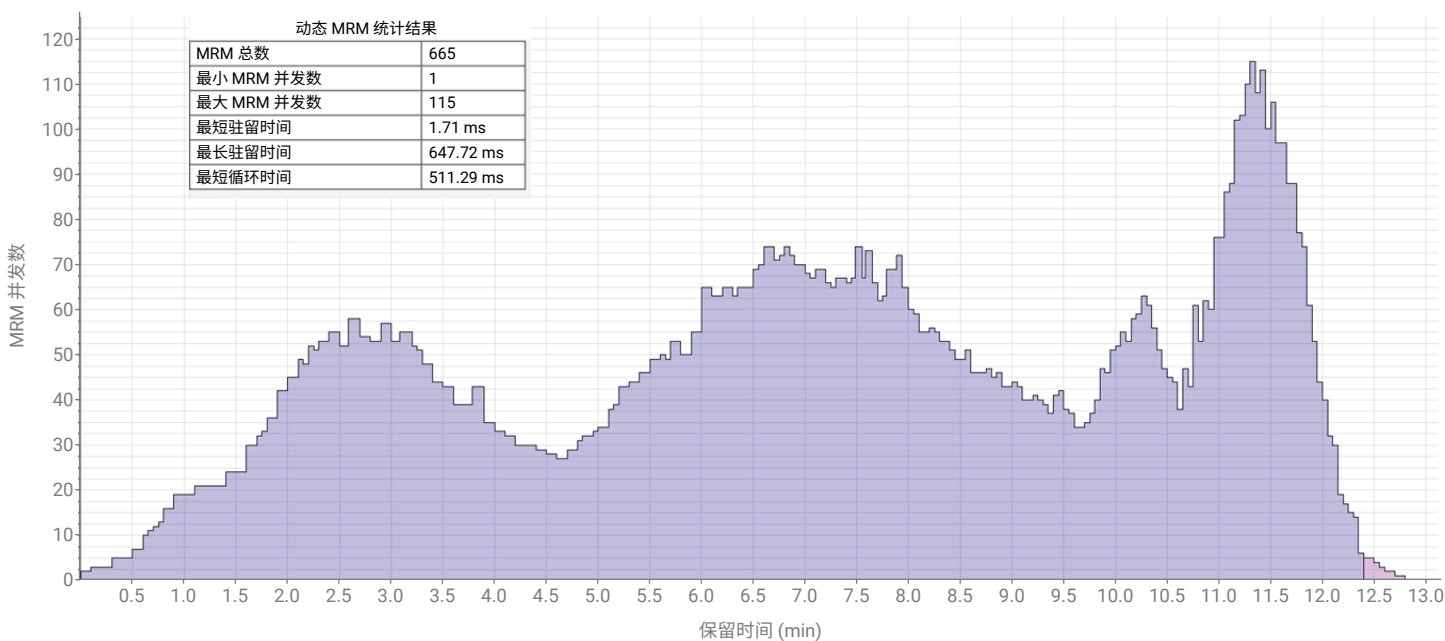


图 3. MRM 并发数与保留时间。显示了各保留时间处的 MRM 并发数的 dMRM 汇总图，以及显示了驻留时间和 MRM 并发数详细信息的 dMRM 统计表



## 血浆中脂质的命名法与结构解析

本文中使用的脂质命名规则遵循 Lipid Maps 联盟制定的指南，以及 Liebisch 等人随后发表的简写符号<sup>[5-7]</sup>。磷脂通常包含两条脂肪酸链，在未进行详细表征的情况下，以碳原子和双键的总组成（即 PC(38:4)）表示。

然而，在确定了酰基链组成的情况下，根据命名规则应将 PC(38:4) 更改为 PC(18:0\_20:4)。这也扩展用于其他脂质类别和亚类。在甘油骨架上脂肪酸的位置已知的情况下，则变为 PC(18:0/20:4)，其中 18:0 在 sn-1 位置，20:4 在 sn-2 位置。已通过色谱分离但尚未完全表征的脂类用 (a) 或 (b) 标记，例如 PC(P-17:0/20:4) (a) 和 PC(P-17:0/20:4) (b)，其中 (a) 和 (b) 表示洗脱顺序。

之前已经报道了该方法基于质量数、碎裂和色谱分析对各种脂质进行结构解析<sup>[1]</sup>。一般情况下，单个甚至多个 MRM 离子对通常无法为给定的脂质种类提供足够的结构信息，尤其是甘油磷脂。有很多方法可以解决这个问题，但每种方法都存在不同的限制。该方法利用离线实验，在特定保留时间提供额外的脂质结构鉴定，并且针对的是具有代表性的混合血浆样品，而不是队列研究中的每个单独样品。参考文献 1 中介绍了此类实验的详细信息，并提供了一些示例：

**磷脂表征：**PC(36:2)，通过离子对 786.6 → 184.1 进行测定，洗脱为两个峰，根据保留时间标记为 PC(36:2) (a) 和 (b)。其他实验使用 PC 类别的正交离子对，在负离子模式下以甲酸盐加合物的形式进行监测，或在正离子模式下使用钠加合物，得到脂肪酸的特定产物。此处，(a) 峰主要为 PC(18:1/18:1)，而第二大的 (b) 峰为 PC(18:0\_18:2)。所有甘油磷脂和鞘磷脂均可以进行此类表征。

**合成标准品：**几种脂质异构体双键的位置有所不同（例如，包含 22:5 脂肪酸的磷脂在血浆中以 ω-3 或 ω-6 异构体的形式存在）。虽然有许多方法可以使用不同的碎裂条件来鉴定这些异构体，但我们选择的是用这两种异构体合成标准品，并通过色谱进行鉴定。同样，可以使用 EID（而不是 CID）等技术或基于洗脱时间鉴定与常规物质（如十七酸）同分异构的单甲基支链脂质（即甲基十六烷酸）。当与血浆样品一起运行时，合成的标准品有助于表征支链和直链酰基链异构体之间的保留时间差异。

**缩醛磷脂：**由于 PC(O) 和 PC(P) 为同分异构体，一种快速确认缩醛磷脂的简单方法是用 HCl 酸雾处理脂质提取物，从而选择性地水解包含烯醚键的物质（缩醛磷脂）。连续运行两个样品，其中一个被水解，即可确认这些物质的身份。

**同位素重叠：**观察到的许多色谱峰可能来自同位素，因为碳以 <sup>13</sup>C 的形式存在，自然丰度约为 1.1%（即，可以在 PC(36:1) 离子对中观察到 PC(36:2) 的 +2 同位素，尽管通常具有不同的保留时间）。这对于磷脂酰胆碱和鞘磷脂尤其成问题，因为它们具有较大的动态范围。通过叠加离子对（在上述示例中为 786.6 → 184.1 和 786.6 → 185.1），可以轻松识别并排除这些物质。如果母离子为同位素，则同位素质量增加只可能发生在子离子中。然后可以排除两个离子对（同位素）中均存在的液相色谱峰。

这些方法共同构成了该方法中结构注释和鉴定的基础。这些实验使用混合血浆样品进行（但也可以使用更常见的参比血浆，例如 NIST1950 血浆）。然后在分析血浆样品时使用这些注释。在使用新基质（例如脑组织）的情况下，由于不同的异构体组成、相对丰度以及血浆中未检测到的新物质/类别，应重复这一过程。此外，本文所介绍的方法侧重于血浆；进一步开发针对其他组织（如大脑）的特定分析将进一步完善此类研究。

## 样品处理和系统配置的注意事项

虽然最终的配置和流程取决于用户，但以下注意事项有助于提高平台的稳定性。

**样品冻融与自动进样器温度：**如果不进行充分控制，脂质在不同温度下的不同溶解度会增加分析的变异性。脂质提取导致血浆脂质被 1:1:0.2 丁醇:甲醇:水混合物所稀释，稀释比为 1:10。在解冻于 -80 °C 下储存的样品时，将其在室温下解冻 1 小时，超声水浴处理 15 分钟，涡旋混合，然后在室温下静置 2 小时后再进样。这可确保脂质在进样前达到平衡。解冻不当会导致脂质浓度随时间发生变化，对于极性较弱的化合物（甘油三酯和胆固醇酯）尤其如此。

**切换阀：**在不损失灵敏度的情况下运行大型样品组的一个关键是，在运行的前 0.8 分钟（含有残留的盐类和极性化合物）切换至废液而不是质谱仪中。这取决于液相色谱的柱体积设置，应根据需要进行调整。

**在线滤芯：**为了延长色谱柱使用寿命并避免堵塞，使用带有 0.3 μm 滤芯（部件号 5067-6189）的在线过滤器可以防止任何沉淀物在系统中积聚。

**单/双柱配置：**通常，如果样品数量较少，单柱配置就已足够，此时每个样品的分析时间约为 16 分钟。但是，如果需要提高通量，双柱配置可以在对一根色谱柱进行再平衡时，用另一根色谱柱进行样品分析，将每个样品的分析时间缩短为 13 分钟。双柱配置的示意图见图 2，两根色谱柱的色谱条件见表 4。

## 峰积分和数据处理

以下是此处详细介绍的实验中所使用的数据处理方案的简要总结。如需了解关于此方法更详细的分步说明，请向安捷伦索取支持文档。

首先将色谱数据导入 Agilent MassHunter 定量分析软件（用于 QQQ）10.0 版或更高版本，并新建一个包含每个待处理样品的批处理文件。然后使用**通过获得的 MRM 数据新建方法**功能生成包含所有相关 MRM 离子对的方法文件，并为标记为内标的每个离子对输入浓度（表 4）。

默认情况下，由 agile2 积分仪进行峰选择，但是，如果积分窗口内包含多个峰，则使用 Spectrum Summation，它允许用户手动选择积分界限。网站 (<https://metabolomics.baker.edu.au/method/lipids>) 中包含所有脂质的注释色谱图，在同分异构体/同质量异构体紧邻洗脱的情况下，可将其作为指南用于确定需要进行积分的峰。注释示例见图 4。如果单个 MRM 离子对有多个需要积分的峰，可以在方法编辑器中复制并重命名该离子对，实现单独积分。



图 4. 网站示例，其中参考了人血浆中要积分的相应峰的内部谱库

将“方法文件”应用到“批处理文件”后，可使用**化合物概览**功能直观地检查所选样品，并确认已选择正确的积分界限。对于保留时间可能因色谱柱性能或溶剂差异而变化的大型批次，由于积分界限固定，Spectrum Summation 将失败。这一问题的示例见图 5。使用时间参比可以解决这

一问题，它根据指定的参比峰自动移动 Spectrum Summation 用于每个 MRM 的定义积分界限（图 5）。如需了解使用时间参比的更多详细说明，以及大型队列的批次比对指南，请向安捷伦索取支持文档。

为了在 MassHunter 定量分析软件中计算浓度（在不使用时间参比的样品组中），需要为每个离子对定义一种内标化合物（表 A2）。应用该方法并完成批处理文件分析后，将自动生成所有样品的最终浓度值。

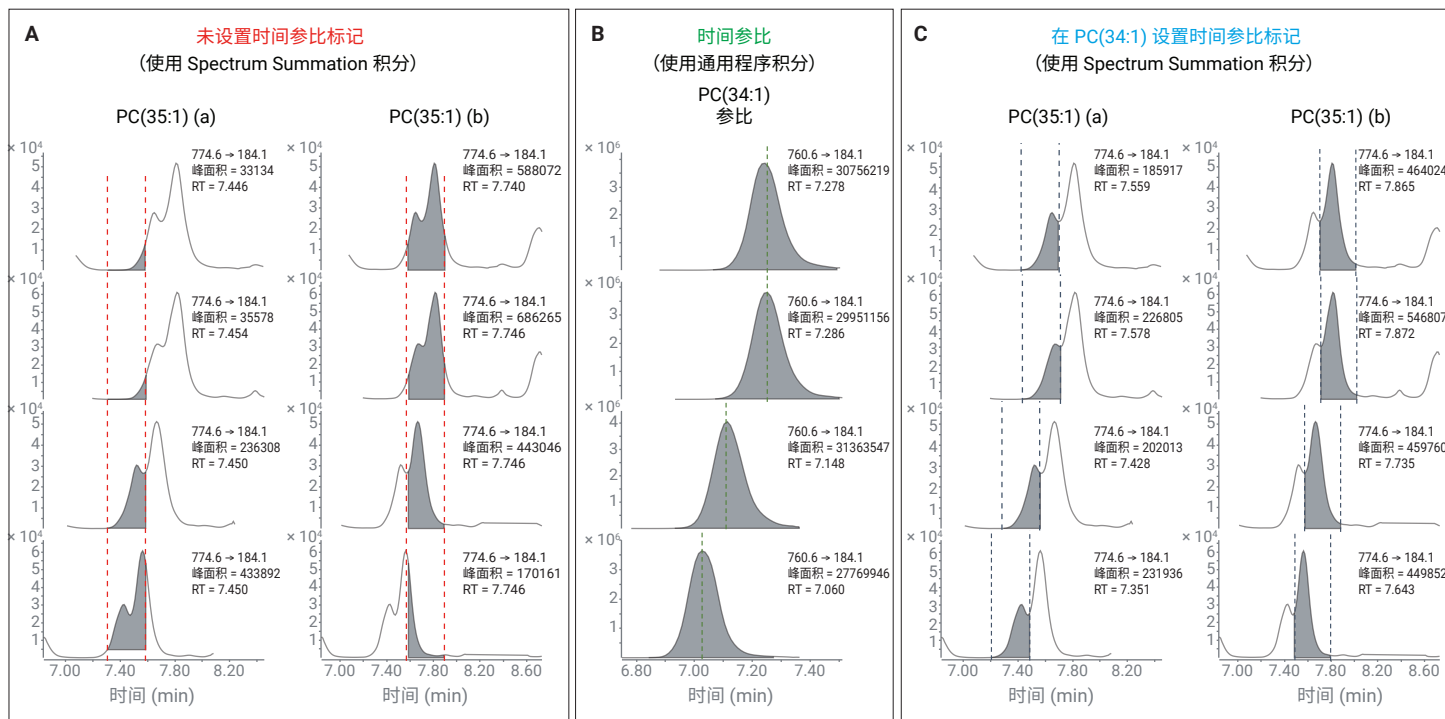


图 5. 使用时间参比化合物比难以积分的峰。(A) PC (35:1) 的两种异构体分别积分，但保留时间漂移导致使用 Spectrum Summation 积分效果不佳。(B) PC(34:1) 丰度高，峰清晰且对称，是一种出色的参比化合物。(C) 启用时间参比并使用 PC(34:1) 作为参比化合物来校正保留时间漂移后，对 PC(35:1) 进行积分

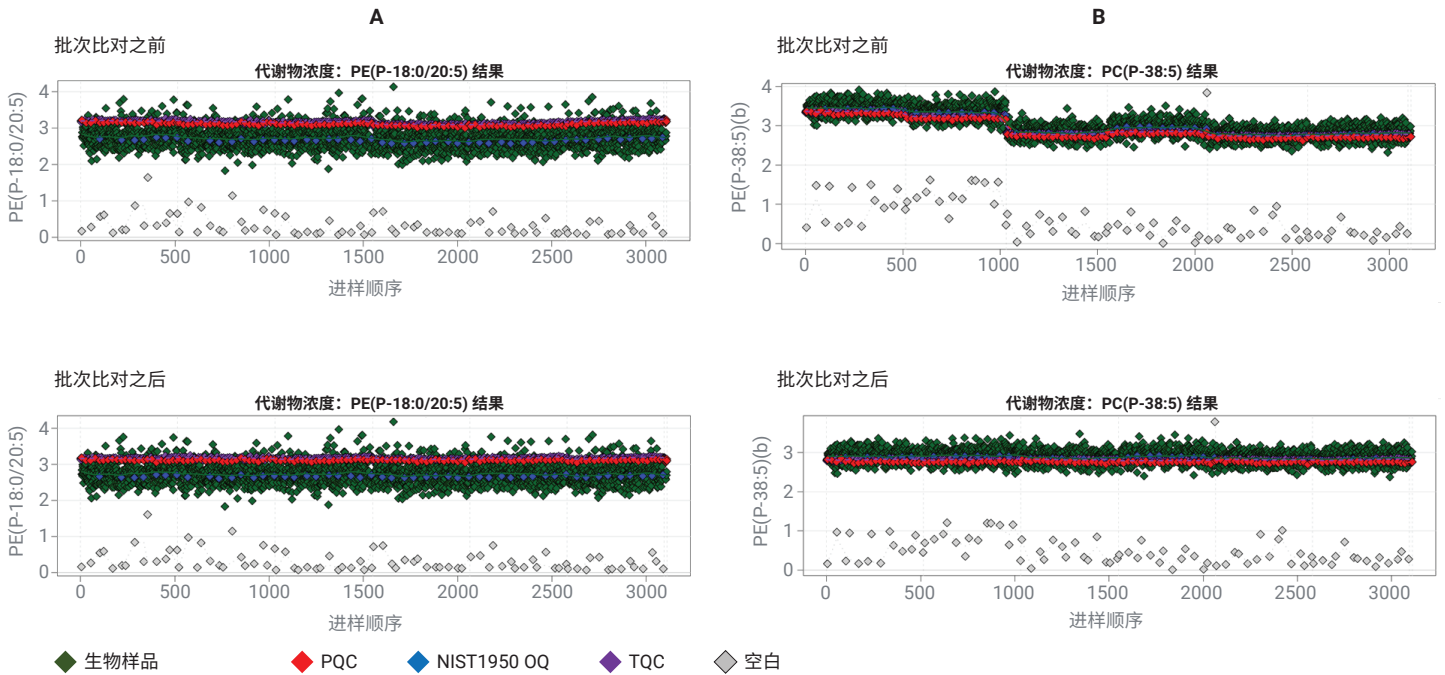


图 6. 使用 PQC 样品对多个批次执行批次比对的示例。(A) 批次之间差异非常小，并且在比对后差异没有增大或改变的脂质示例。(B) 不同批次之间差异较大的脂质示例，以及比对后的改善。所有图中均显示 6 个批次共 2500 个生物样品 (X 轴)，以及用 log<sub>2</sub> 对数值表示的脂质丰度 (Y 轴)

## 结果与讨论

为了确定该方法的重现性，依次进样 50 个 TQCs (参见图 7 的色谱图示例) 并计算了峰面积和脂质浓度的 %CV。763 种脂质的原始峰面积的 %CV 平均值和中值分别为 5.1% 和 4.1%。计算每种脂质的浓度后，%CV 的平均值和中值分别为 5.6% 和 4.0%。每种脂质各自的 %CV 见表 A2。对于这些脂质，应考虑通常由萃取和液相色谱分析过程引入的溶剂和增塑剂产生的背景信号。表 A2 显示了每种脂质的背景水平，以 NIST1950 QC 样品产生的信号的百分比表示。585 种脂质的背景 < 1.0%，仅 28 种脂质的背景 > 20.0%。应特别注意包含饱和脂肪酸链的游离脂肪酸和二酰甘油，以及一些丰度极低的脂质种类。

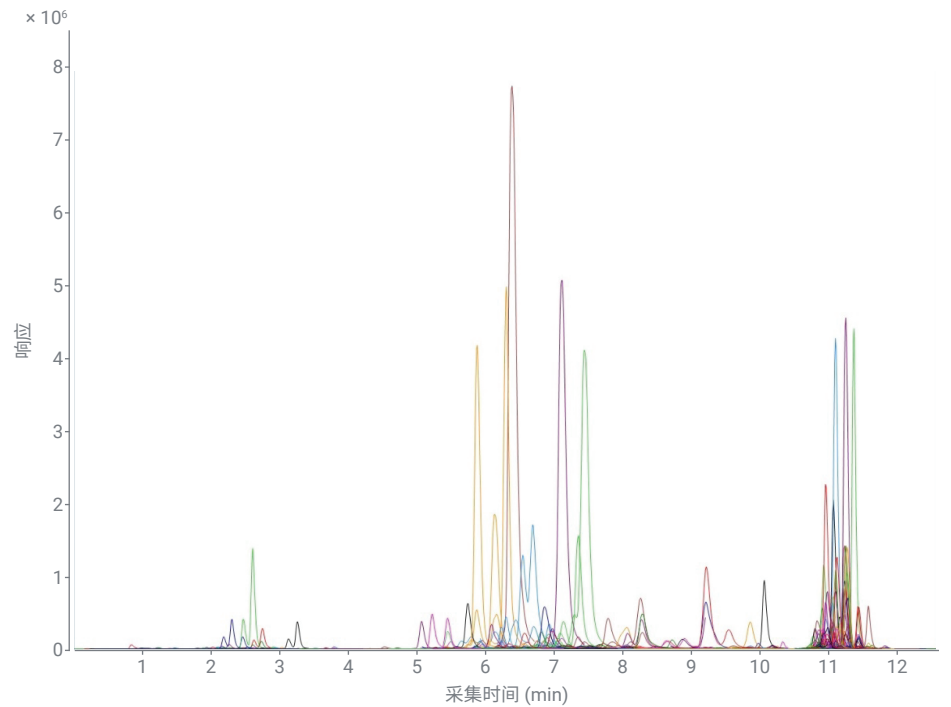


图 7. 用此方法测得的脂质的叠加 MRM 色谱图

通过运行包含 2500 个血浆样品的队列，进一步证明了方法的稳定性。该运行分为六个批次，每个批次包含 417 个样品、9 个 NIST1950 QCs、23 个 PQCs、14 个空白和 25 个 TQCs。单个代表性批次和所有比对批次的脂质浓度 %CV 汇总见表 7。对于每个质控样品，图 8 显示了 CV% 直方图，突出显示了处于 CV 阈值范围内的脂质数量。在质控样品中，约有 600 种脂质的 CV 小于 10%。然后对所有 QCs 和样品进行主成分分析（图 9），以突出显示脂质相对于质控样品的生物学差异。

本应用简报详细介绍了一种稳定且高度优化的 LC/MS/MS 方案，用于测定血浆样品或同类物质中的 750 多种不同脂质。该方案运行时间短、数据重现性高，并且能够在多个批次甚至不同实验室之间进行数据比对，有助于整合大型脂质组数据集。这些数据对于阐明脂质种类与疾病转归之间的关联，从而更好地了解疾病发病机制至关重要。

表 7. 单个批次以及多个批次中每种 QC 类型的 %CV

	NIST 1950 QC		PQC		TQC	
	均值	中值	均值	中值	均值	中值
单个批次 (417 个样品)	9.7	7.7	7.9	6.6	7.0	5.6
所有批次 (2500 个样品)	9.7	7.7	8.9	7.4	8.6	7.0

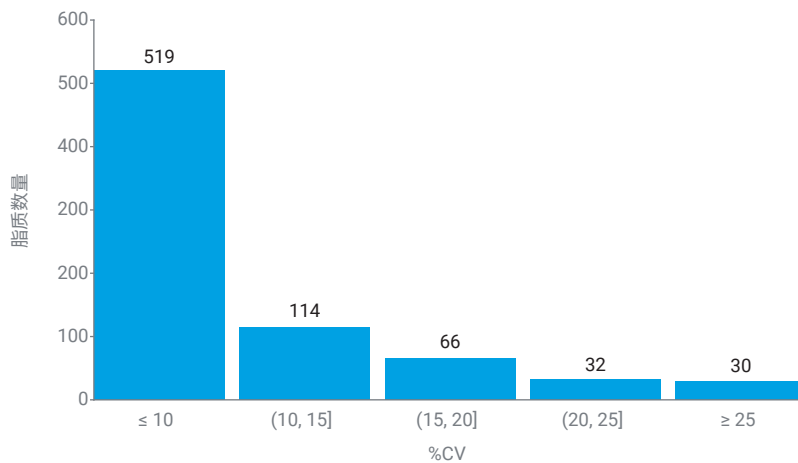


图 8. 在一个批次中测量的各种 QC 样品中脂质变异系数的直方图

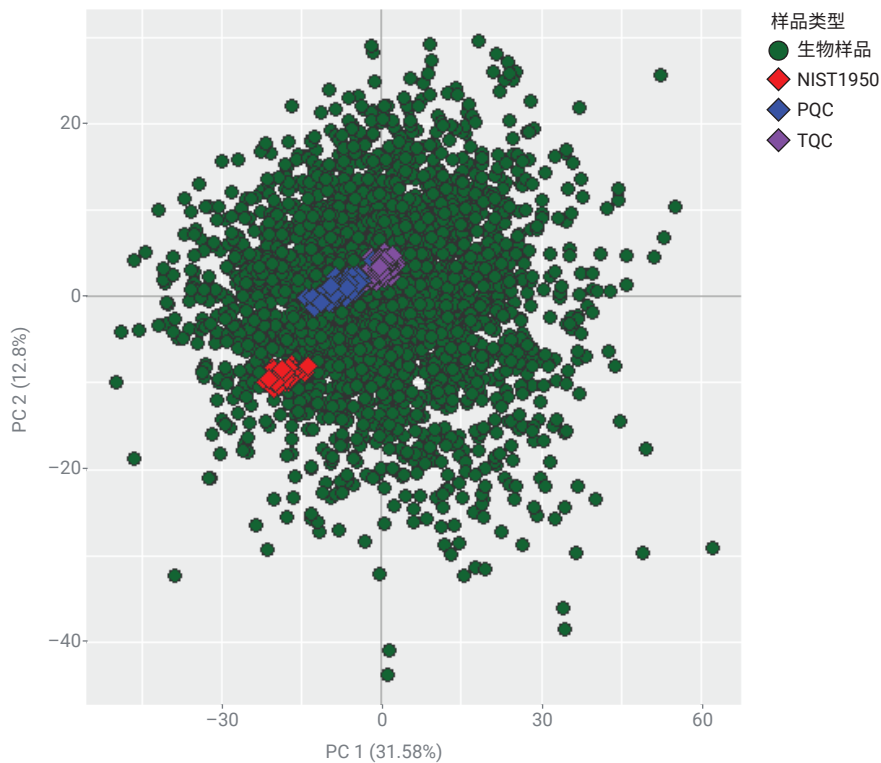


图 9. 来自一批 2500 个生物样品和所有相关 QC 样品的脂质组的 PCA 结果

该方法的一个局限性在于一些异构体不能完全分离，而这必然受到了运行时间有限的限制。但请注意，通过正交实验、使用各种加合物或电离模式，或扩展色谱运行可以突破这些限制。对应地，用户必须在分析时间和特异性方面做出妥协。

LC/MS/MS 采集脂质组学定量数据的能力在领域内仍然是一个存在争议的话题。事实上，许多因素会对数据的定量准确度造成影响，其中最为显著的是基质效应、差异抑制以及各种脂质的不同离子化效率。对于绝对定量，还必须校正标准品和目标分析物之间的同位素分布差异。出于这些原因，该方法被称为相对定量，所获得的值在样品之间具有可比性，但不一定是样品的绝对浓度。可以合理地认为，在大多数情况下，研究目的在于确定关联性（即随年龄变化的脂质代谢物）或样品组（健康组与糖尿病组）之间的倍数差异。在这类分析中，绝对定量与此处获得的结果相比没有任何附加价值。当需要在研究之间进行比较时（绝对定量具有一定的优势），使用本文所述的标准物质（例如 NIST SRM 1950）进行比对可以解决这个问题。

该方案涵盖广泛的脂质代谢物，并且不可能为每种化合物都提供一种内标。因此，大多数脂质类别使用同一个内标。尽管存在这些限制，该方法在每种脂质化合物的相对定量中仍然具有良好的重现性和精密

度。事实上，大多数研究的目标是确定样品之间脂质浓度的相对差异，或确定人群中脂质种类与参与者特征或临床结果之间的关联性。在这种情况下，脂质的绝对定量除了用于比较不同平台之间的结果外，几乎没有任何作用。

在需要对不同研究或实验室之间的数据进行比较的情况下，安捷伦会使用所有实验中通常包含的 NIST SRM 1950 血浆 QC 样品。该样品可以从多个供应商处轻松获得，并为所有测量的脂质提供了一个共同的参考点。因此，只要脂质测量处于仪器的线性动态范围内，就可以使用 NIST QC 样品进行标准化和比较。

该方案可轻松调整以适应其他生物样品。然而，考虑到组织脂质组之间的潜在差异，建议使用上述方法对新的组织类型进行组织特异性注释。在未进行组织特异性注释的情况下，建议在检测其他基质时将注释恢复到总组成水平（例如，PC(36:2) (a) 和 (b)）。组织特异性注释可能包含组织特异性脂质，例如心磷脂，它是一种仅存在于线粒体中的脂质，在血浆或所述方法中不存在。还可以对色谱进行其他调整，尤其是在不要求高通量的小型研究中，因为更长、更平缓的梯度能够进一步改善脂质分离（包括同分异构体和同质量异构体），从而分离更多的脂质异构体，减少来自共洗脱物质的信号抑制，并进一步改善每个离子对的驻留时间。

## 参考文献

1. Huynh, K. *et al.* High-Throughput Plasma Lipidomics: Detailed Mapping of the Associations with Cardiometabolic Risk Factors. *Cell Chem. Biol.* **2019**, *26*(1), 71–84.e4
2. Beyene, H. B. *et al.* High-Coverage Plasma Lipidomics Reveals Novel Sex-Specific Lipidomic Fingerprints of Age and BMI: Evidence from Two Large Population Cohort Studies. *PLoS Biol.* **2020**, *18*(9), e3000870
3. Alshehry, Z. H. *et al.* An Efficient Single Phase Method for the Extraction of Plasma Lipids. *Metabolites* **2015**, *5*(2), 389–403
4. Muniandy, M. 等，采用 Agilent Bravo 自动化液体处理平台的半自动化脂质萃取方案，*安捷伦科技公司应用简报*，出版号 5991-5724CHCN，**2015**
5. Liebisch, G. *et al.* Shorthand Notation for Lipid Structures Derived from Mass Spectrometry. *J. Lipid Res.* **2013**, *54*(6), 1523–1530
6. Fahy, E. *et al.* A Comprehensive Classification System for Lipids. *J. Lipid Res.* **2005**, *46*(5), 839–861
7. Fahy, E. *et al.* Update of the LIPID MAPS Comprehensive Classification System for Lipids. *J. Lipid Res.* **2009**, *50*, S9–14

## 附录

表 A1. 所有 dMRM 离子对的 Agilent MassHunter 采集软件详细信息<sup>1</sup>

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1 分辨率	子离子	MS2 分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
1	AcylCarn	AC(10:0)	FALSE	316.3	Unit	85.1	Unit	0.9	0.6	166	30	5	Positive
2	AcylCarn	AC(12:0)	FALSE	344.3	Unit	85.1	Unit	1.2	0.6	166	30	5	Positive
3	AcylCarn	AC(12:1)	FALSE	342.3	Unit	85.1	Unit	1.0	0.6	166	30	5	Positive
4	AcylCarn	AC(13:0)	FALSE	358.3	Unit	85.1	Unit	1.4	0.6	166	30	5	Positive
5	AcylCarn	AC(14:0)	FALSE	372.3	Unit	85.1	Unit	1.8	0.6	166	30	5	Positive
6	AcylCarn	AC(14:0)-OH	FALSE	388.3	Unit	85.1	Unit	1.4	0.6	166	30	5	Positive
7	AcylCarn	AC(14:1)	FALSE	370.3	Unit	85.1	Unit	1.5	0.6	166	30	5	Positive
8	AcylCarn	AC(14:1)-OH	FALSE	386.3	Unit	85.1	Unit	1.1	0.6	166	30	5	Positive
9	AcylCarn	AC(14:2)	FALSE	368.3	Unit	85.1	Unit	1.2	0.6	166	30	5	Positive
10	AcylCarn	AC(15:0) (a\b)	FALSE	386.3	Unit	85.1	Unit	2.1	1	166	30	5	Positive
11	AcylCarn	AC(16:0)	FALSE	400.4	Unit	85.1	Unit	2.5	0.6	166	30	5	Positive
12	AcylCarn	AC(16:0)-OH	FALSE	416.4	Unit	85.1	Unit	2.0	0.6	166	30	5	Positive
13	AcylCarn	AcylCarnitine 16:0 d3 (IS)	TRUE	403.4	Unit	85.1	Unit	2.5	0.6	166	30	5	Positive
14	AcylCarn	AC(16:1)	FALSE	398.3	Unit	85.1	Unit	2.0	0.6	166	30	5	Positive
15	AcylCarn	AC(16:1)-OH	FALSE	414.3	Unit	85.1	Unit	1.6	0.6	166	30	5	Positive
16	AcylCarn	AC(17:0) (a\b)	FALSE	414.4	Unit	85.1	Unit	2.8	0.6	166	30	5	Positive
17	AcylCarn	AC(18:0)	FALSE	428.4	Unit	85.1	Unit	3.2	0.6	166	30	5	Positive
18	AcylCarn	AC(18:0)-OH	FALSE	444.4	Unit	85.1	Unit	2.7	0.6	166	30	5	Positive
19	AcylCarn	AC(18:1)	FALSE	426.4	Unit	85.1	Unit	2.7	0.6	166	30	5	Positive
20	AcylCarn	AC(18:1)-OH	FALSE	442.4	Unit	85.1	Unit	2.2	0.6	166	30	5	Positive
21	AcylCarn	AC(18:2)	FALSE	424.3	Unit	85.1	Unit	2.2	0.6	166	30	5	Positive
22	AcylCarn	AC(18:3)	FALSE	422.3	Unit	85.1	Unit	1.9	0.6	166	30	5	Positive
23	AcylCarn	AC(20:3) (a\b)	FALSE	450.3	Unit	85.1	Unit	2.5	0.6	166	30	5	Positive
24	AcylCarn	AC(20:3)-OH	FALSE	466.3	Unit	85.1	Unit	2.1	0.6	166	30	5	Positive
25	AcylCarn	AC(20:4)	FALSE	448.3	Unit	85.1	Unit	2.2	0.6	166	30	5	Positive
26	AcylCarn	AC(20:5)	FALSE	446.3	Unit	85.1	Unit	1.8	0.6	166	30	5	Positive
27	AcylCarn	AC(22:5)	FALSE	474.3	Unit	85.1	Unit	2.3	0.6	166	30	5	Positive
28	AcylCarn	AC(22:5)-OH	FALSE	490.3	Unit	85.1	Unit	2.0	0.6	166	30	5	Positive
29	AcylCarn	AC(22:6)	FALSE	472.3	Unit	85.1	Unit	2.2	0.6	166	30	5	Positive
30	AcylCarn	AC(24:0)	FALSE	512.3	Unit	85.1	Unit	5.0	0.6	166	30	5	Positive
31	AcylCarn	AC(24:1) (a\b)	FALSE	510.3	Unit	85.1	Unit	4.4	0.6	166	30	5	Positive
32	AcylCarn	AC(26:0)	FALSE	540.3	Unit	85.1	Unit	5.9	0.6	166	30	5	Positive
33	AcylCarn	AC(26:1)	FALSE	538.3	Unit	85.1	Unit	5.0	0.6	166	30	5	Positive
34	CE	CE(14:0)	FALSE	614.6	Unit	369.3	Unit	11.6	0.4	166	10	5	Positive
35	CE	CE(15:0)	FALSE	628.6	Unit	369.3	Unit	11.7	0.4	166	10	5	Positive
36	CE	CE(16:0)	FALSE	642.6	Unit	369.3	Unit	11.8	0.4	166	10	5	Positive
37	CE	CE(16:1)	FALSE	640.6	Unit	369.3	Unit	11.6	0.4	166	10	5	Positive
38	CE	CE(16:2)	FALSE	638.6	Unit	369.3	Unit	11.4	0.4	166	10	5	Positive
39	CE	CE(17:0)	FALSE	656.6	Unit	369.3	Unit	11.9	0.4	166	10	5	Positive
40	CE	CE(17:1)	FALSE	654.6	Unit	369.3	Unit	11.7	0.4	166	10	5	Positive
41	CE	CE(18:0)	FALSE	670.7	Unit	369.3	Unit	12.1	0.4	166	10	5	Positive
42	CE	CE 18:0-d6 (IS)	TRUE	676.7	Unit	375.3	Unit	12.0	0.4	166	10	5	Positive
43	CE	CE(18:1)	FALSE	668.6	Unit	369.3	Unit	11.8	0.4	166	10	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
44	CE	CE(18:2)	FALSE	666.6	Unit	369.3	Unit	11.6	0.4	166	10	5	Positive
45	oxLipid	CE(18:2) [+OH]	FALSE	682.6	Unit	369.3	Unit	10.9	0.6	166	10	5	Positive
46	CE	CE(18:3)	FALSE	664.6	Unit	369.3	Unit	11.4	0.4	166	10	5	Positive
47	CE	CE(20:0)	FALSE	698.7	Unit	369.3	Unit	12.2	0.6	166	10	5	Positive
48	CE	CE(20:1)	FALSE	696.7	Unit	369.3	Unit	12.0	0.4	166	10	5	Positive
49	CE	CE(20:2)	FALSE	694.7	Unit	369.3	Unit	11.8	0.4	166	10	5	Positive
50	CE	CE(20:3)	FALSE	692.6	Unit	369.3	Unit	11.7	0.4	166	10	5	Positive
51	CE	CE(20:4)	FALSE	690.6	Unit	369.3	Unit	11.5	0.4	166	10	5	Positive
52	oxLipid	CE(20:4) [+OH]	FALSE	706.6	Unit	369.3	Unit	10.9	0.6	166	10	5	Positive
53	CE	CE(20:5)	FALSE	688.6	Unit	369.3	Unit	11.3	0.4	166	10	5	Positive
54	CE	CE(22:0)	FALSE	726.7	Unit	369.3	Unit	12.4	0.6	166	10	5	Positive
55	CE	CE(22:1)	FALSE	724.7	Unit	369.3	Unit	12.2	0.4	166	10	5	Positive
56	CE	CE(22:4)	FALSE	718.7	Unit	369.3	Unit	11.7	0.4	166	10	5	Positive
57	CE	CE(22:5)	FALSE	716.6	Unit	369.3	Unit	11.5	1	166	10	5	Positive
58	CE	CE(22:6)	FALSE	714.6	Unit	369.3	Unit	11.4	0.4	166	10	5	Positive
59	CE	CE(24:0)	FALSE	754.7	Unit	369.3	Unit	12.6	0.4	166	10	5	Positive
60	CE	CE(24:1)	FALSE	752.7	Unit	369.3	Unit	12.4	0.4	166	10	5	Positive
61	CE	CE(24:4)	FALSE	746.7	Unit	369.3	Unit	11.9	0.4	166	10	5	Positive
62	CE	CE(24:5)	FALSE	744.7	Unit	369.3	Unit	11.7	0.4	166	10	5	Positive
63	CE	CE(24:6)	FALSE	742.7	Unit	369.3	Unit	11.5	0.4	166	10	5	Positive
64	Cer1P	Cer1P(d18:1/16:0)	FALSE	618.424	Unit	264.3	Unit	5.5	0.6	166	29	4	Positive
65	Cer	Cer(d16:1/16:0)	FALSE	510.6	Unit	236.3	Unit	6.7	0.5	166	25	5	Positive
66	Cer	Cer(d16:1/18:0)	FALSE	538.6	Unit	236.3	Unit	7.8	0.5	166	25	5	Positive
67	Cer	Cer(d16:1/20:0)	FALSE	566.6	Unit	236.3	Unit	9.1	0.5	166	25	5	Positive
68	Cer	Cer(d16:1/22:0)	FALSE	594.6	Unit	236.3	Unit	10.1	0.5	166	25	5	Positive
69	Cer	Cer(d16:1/23:0)	FALSE	608.6	Unit	236.3	Unit	10.2	0.5	166	25	5	Positive
70	Cer	Cer(d16:1/24:0)	FALSE	622.6	Unit	236.3	Unit	10.3	0.5	166	25	5	Positive
71	Cer	Cer(d16:1/24:1)	FALSE	620.6	Unit	236.3	Unit	10.1	0.5	166	25	5	Positive
72	Cer	Cer(d17:1/16:0)	FALSE	524.6	Unit	250.3	Unit	7.2	0.5	166	25	5	Positive
73	Cer	Cer(d17:1/18:0)	FALSE	552.6	Unit	250.3	Unit	8.4	0.5	166	25	5	Positive
74	Cer	Cer(d17:1/20:0)	FALSE	580.6	Unit	250.3	Unit	9.7	0.5	166	25	5	Positive
75	Cer	Cer(d17:1/22:0)	FALSE	608.6	Unit	250.3	Unit	10.2	0.5	166	25	5	Positive
76	Cer	Cer(d17:1/23:0)	FALSE	622.6	Unit	250.3	Unit	10.3	0.5	166	25	5	Positive
77	Cer	Cer(d17:1/24:0)	FALSE	636.6	Unit	250.3	Unit	10.4	0.5	166	25	5	Positive
78	Cer	Cer(d17:1/24:1)	FALSE	634.6	Unit	250.3	Unit	10.2	0.5	166	25	5	Positive
79	Cer	Cer(d18:1/16:0)	FALSE	538.5	Unit	264.3	Unit	7.8	0.7	166	25	5	Positive
80	Cer	Cer(d18:1/18:0)	FALSE	566.6	Unit	264.3	Unit	9.0	0.7	166	25	5	Positive
81	Cer	Cer(d18:1/20:0)	FALSE	594.6	Unit	264.3	Unit	10.1	0.7	166	25	5	Positive
82	Cer	Cer(d18:1/21:0)	FALSE	608.6	Unit	264.3	Unit	10.2	0.7	166	25	5	Positive
83	Cer	Cer(d18:1/22:0)	FALSE	622.6	Unit	264.3	Unit	10.3	0.7	166	25	5	Positive
84	Cer	Cer(d18:1/23:0)	FALSE	636.6	Unit	264.3	Unit	10.4	0.7	166	25	5	Positive
85	Cer	Cer(d18:1/24:0)	FALSE	650.6	Unit	264.3	Unit	10.5	0.7	166	25	5	Positive
86	Cer	Cer(d18:1/24:1)	FALSE	648.6	Unit	264.3	Unit	10.3	0.7	166	25	5	Positive
87	Cer	Cer(d18:1/26:0)	FALSE	678.6	Unit	264.3	Unit	10.7	0.7	166	25	5	Positive
88	Cer	Cer(d18:1-d7/18:0) (IS)	TRUE	573.6	Unit	271.4	Unit	9.0	1	166	25	5	Positive
89	Cer	Cer(d18:2/16:0)	FALSE	536.5	Unit	262.3	Unit	6.9	0.7	166	25	5	Positive
90	Cer	Cer(d18:2/18:0)	FALSE	564.6	Unit	262.3	Unit	8.1	0.7	166	25	5	Positive



化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
91	Cer	Cer(d18:2/20:0)	FALSE	592.6	Unit	262.3	Unit	9.3	0.5	166	25	5	Positive
92	Cer	Cer(d18:2/22:0)	FALSE	620.6	Unit	262.3	Unit	10.1	0.5	166	25	5	Positive
93	Cer	Cer(d18:2/23:0)	FALSE	634.6	Unit	262.3	Unit	10.2	0.5	166	25	5	Positive
94	Cer	Cer(d18:2/24:0)	FALSE	648.6	Unit	262.3	Unit	10.4	0.5	166	25	5	Positive
95	Cer	Cer(d18:2/24:1)	FALSE	646.6	Unit	262.3	Unit	10.1	0.5	166	25	5	Positive
96	Cer	Cer(d18:2/26:0)	FALSE	676.6	Unit	262.3	Unit	10.5	0.5	166	25	5	Positive
97	Cer	Cer(d19:1/18:0)	FALSE	580.6	Unit	278.3	Unit	9.4	0.5	166	25	5	Positive
98	Cer	Cer(d19:1/20:0)	FALSE	608.6	Unit	278.3	Unit	10.1	0.5	166	25	5	Positive
99	Cer	Cer(d19:1/22:0)	FALSE	636.6	Unit	278.3	Unit	10.4	0.5	166	25	5	Positive
100	Cer	Cer(d19:1/23:0)	FALSE	650.6	Unit	278.3	Unit	10.5	0.5	166	25	5	Positive
101	Cer	Cer(d19:1/24:0)	FALSE	664.6	Unit	278.3	Unit	10.5	0.5	166	25	5	Positive
102	Cer	Cer(d19:1/24:1)	FALSE	662.6	Unit	278.3	Unit	10.4	0.5	166	25	5	Positive
103	Cer	Cer(d20:1/22:0)	FALSE	650.6	Unit	292.3	Unit	10.5	0.5	166	25	5	Positive
104	Cer	Cer(d20:1/23:0)	FALSE	664.6	Unit	292.3	Unit	10.6	0.5	166	25	5	Positive
105	Cer	Cer(d20:1/24:0)	FALSE	678.6	Unit	292.3	Unit	10.7	0.5	166	25	5	Positive
106	Cer	Cer(d20:1/24:1)	FALSE	676.6	Unit	292.3	Unit	10.5	0.5	166	25	5	Positive
107	m18:0	Cer(m18:0/20:0)	FALSE	580.6	Unit	268.4	Unit	10.3	0.6	166	35	5	Positive
108	m18:0	Cer(m18:0/22:0)	FALSE	608.6	Unit	268.4	Unit	10.5	0.6	166	35	5	Positive
109	m18:0	Cer(m18:0/23:0)	FALSE	622.6	Unit	268.4	Unit	10.6	0.6	166	35	5	Positive
110	m18:0	Cer(m18:0/24:0)	FALSE	636.6	Unit	268.4	Unit	10.7	0.6	166	35	5	Positive
111	m18:0	Cer(m18:0/24:1)	FALSE	634.6	Unit	268.4	Unit	10.5	0.6	166	35	5	Positive
112	m18:1	Cer(m18:1/18:0)	FALSE	550.6	Unit	266.4	Unit	9.3	0.6	166	35	5	Positive
113	m18:1	Cer(m18:1/20:0)	FALSE	578.6	Unit	266.4	Unit	10.1	0.6	166	35	5	Positive
114	m18:1	Cer(m18:1/22:0)	FALSE	606.6	Unit	266.4	Unit	10.3	0.6	166	35	5	Positive
115	m18:1	Cer(m18:1/23:0)	FALSE	620.6	Unit	266.4	Unit	10.4	0.6	166	35	5	Positive
116	m18:1	Cer(m18:1/24:0)	FALSE	634.6	Unit	266.4	Unit	10.5	0.6	166	35	5	Positive
117	m18:1	Cer(m18:1/24:1)	FALSE	632.6	Unit	266.4	Unit	10.4	0.6	166	35	5	Positive
118	Bile Acids	CA	FALSE	426.3	Unit	355.3	Unit	1.0	2	166	19	4	Positive
119	Bile Acids	Cholic Acid d4 (IS)	TRUE	430.3	Unit	359.3	Unit	1.0	1	166	19	4	Positive
120	COH	COH	FALSE	369.4	Unit	161.2	Unit	6.4	0.6	166	19	5	Positive
121	COH	COH-d7 (IS)	TRUE	376.4	Unit	161.2	Unit	6.4	0.6	166	19	5	Positive
122	DE	DE(16:0)	FALSE	640.6	Unit	367.4	Unit	11.6	0.6	166	10	4	Positive
123	DE	DE(18:1)	FALSE	666.6	Unit	367.4	Unit	11.6	0.6	166	10	4	Positive
124	DE	DE(18:1) ester d6 (IS)	TRUE	672.6	Unit	373.4	Unit	11.6	0.6	166	10	4	Positive
125	DE	DE(18:2)	FALSE	664.6	Unit	367.4	Unit	11.4	0.6	166	10	4	Positive
126	DE	DE(20:4)	FALSE	688.6	Unit	367.4	Unit	11.3	0.6	166	10	4	Positive
127	Bile Acids	dxCA	FALSE	410.3	Unit	357.3	Unit	1.6	2	166	15	5	Positive
128	DG	DG(15:0 18:1) d7 (IS)	TRUE	605.5	Unit	299.5	Unit	9.6	1	166	25	4	Positive
129	DG	DG(14:0_16:0)	FALSE	558.5	Unit	313.3	Unit	8.8	1.4	166	25	5	Positive
130	DG	DG(16:0_16:0)	FALSE	586.5	Unit	313.2	Unit	10.1	1.4	166	25	5	Positive
131	DG	DG(16:0_16:1)	FALSE	584.5	Unit	313.2	Unit	9.1	1.4	166	25	5	Positive
132	DG	DG(14:0_18:2)	FALSE	582.5	Unit	285.2	Unit	8.1	1.4	166	25	5	Positive
133	DG	DG(16:0_18:1)	FALSE	612.6	Unit	313.3	Unit	10.0	1.4	166	25	5	Positive
134	DG	DG(16:1_18:1)	FALSE	610.5	Unit	339.2	Unit	9.3	1.4	166	25	5	Positive
135	DG	DG(16:0_18:2)	FALSE	610.5	Unit	313.2	Unit	9.4	1.4	166	25	5	Positive
136	DG	DG(18:0_18:1)	FALSE	640.6	Unit	341.3	Unit	10.2	1.4	166	25	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
137	DG	DG(18:1_18:1)	FALSE	638.6	Unit	339.3	Unit	10.1	1.4	166	25	5	Positive
138	DG	DG(18:0_18:2)	FALSE	638.6	Unit	341.3	Unit	10.1	1.4	166	25	5	Positive
139	DG	DG(18:1_18:2)	FALSE	636.6	Unit	339.3	Unit	9.6	1.4	166	25	5	Positive
140	DG	DG(18:2_18:2)	FALSE	634.5	Unit	337.2	Unit	8.7	1.4	166	25	5	Positive
141	DG	DG(18:1_18:3)	FALSE	634.5	Unit	339.2	Unit	8.8	1.4	166	25	5	Positive
142	DG	DG(16:0_20:4)	FALSE	634.5	Unit	313.2	Unit	9.2	1.4	166	25	5	Positive
143	DG	DG(18:1_20:3)	FALSE	662.6	Unit	339.3	Unit	10.0	1.4	166	25	5	Positive
144	DG	DG(18:0_20:4)	FALSE	662.6	Unit	341.3	Unit	10.1	1.4	166	25	5	Positive
145	DG	DG(18:1_20:4)	FALSE	660.6	Unit	339.3	Unit	9.4	1.4	166	25	5	Positive
146	DG	DG(16:0_22:5)	FALSE	660.6	Unit	313.3	Unit	9.3	1.4	166	25	5	Positive
147	DG	DG(18:2_20:4)	FALSE	658.5	Unit	337.2	Unit	8.5	1.4	166	25	5	Positive
148	DG	DG(16:0_22:6)	FALSE	658.5	Unit	313.2	Unit	8.9	1.4	166	25	5	Positive
149	DG	DG(18:1_20:5)	FALSE	658.6	Unit	339.3	Unit	8.6	1.4	166	25	5	Positive
150	DG	DG(18:1_22:5)	FALSE	686.6	Unit	339.3	Unit	9.8	1.6	166	25	5	Positive
151	DG	DG(18:1_22:6)	FALSE	684.6	Unit	339.3	Unit	9.1	1.4	166	25	5	Positive
152	DG	DG(18:2_22:6)	FALSE	682.6	Unit	337.3	Unit	8.2	1.4	166	25	5	Positive
153	dhCer	dhCer(d18:0/16:0)	FALSE	540.5	Unit	284.3	Unit	8.3	0.6	166	27	4	Positive
154	dhCer	dhCer(d18:0/18:0)	FALSE	568.6	Unit	284.3	Unit	9.8	0.8	166	27	4	Positive
155	dhCer	dhCer(d18:0/20:0)	FALSE	596.6	Unit	284.3	Unit	10.2	0.8	166	27	4	Positive
156	dhCer	dhCer(d18:0/22:0)	FALSE	624.6	Unit	284.3	Unit	10.4	0.8	166	27	4	Positive
157	dhCer	dhCer(d18:0/24:0)	FALSE	652.7	Unit	284.3	Unit	10.6	0.8	166	27	4	Positive
158	dhCer	dhCer(d18:0/24:1)	FALSE	650.6	Unit	284.3	Unit	10.4	0.8	166	27	4	Positive
159	dhCer	dhCer 8:0 (IS)	TRUE	428.4	Unit	284.3	Unit	4.6	0.8	166	27	4	Positive
160	dimethyl-CE	dimethyl-CE(18:1)	FALSE	696.6	Unit	397.3	Unit	12.0	0.4	166	10	4	Positive
161	dimethyl-CE	dimethyl-CE(18:2)	FALSE	694.6	Unit	397.3	Unit	11.8	0.4	166	10	4	Positive
162	dimethyl-CE	dimethyl-CE(20:4)	FALSE	718.6	Unit	397.3	Unit	11.7	0.4	166	10	4	Positive
163	FFA	FA(14:0)	FALSE	227.2	Unit	227.2	Unit	3.1	2	166	0	4	Negative
164	FFA	FA(16:0)	FALSE	255.2	Unit	255.2	Unit	3.8	2	166	0	4	Negative
165	FFA	FA(16:1)	FALSE	253.2	Unit	253.2	Unit	3.3	2	166	0	4	Negative
166	FFA	FA(17:0)	FALSE	269.3	Unit	269.2	Unit	4.1	2	166	0	4	Negative
167	FFA	FA(17:1)	FALSE	267.2	Unit	267.2	Unit	3.6	2	166	0	4	Negative
168	FFA	FA(18:0)	FALSE	283.3	Unit	283.3	Unit	4.4	2	166	0	4	Negative
169	FFA	FA(18:1)	FALSE	281.3	Unit	281.2	Unit	3.9	2	166	0	4	Negative
170	FFA	FA(18:1) d9 (IS)	TRUE	290.3	Unit	290.2	Unit	3.9	2	166	0	4	Negative
171	FFA	FA(18:2)	FALSE	279.2	Unit	279.2	Unit	3.5	2	166	0	4	Negative
172	FFA	FA(18:3)	FALSE	277.2	Unit	277.2	Unit	3.1	2	166	0	4	Negative
173	FFA	FA(20:2)	FALSE	307.3	Unit	307.3	Unit	4.1	2	166	0	4	Negative
174	FFA	FA(20:3)	FALSE	305.3	Unit	305.2	Unit	3.8	2	166	0	4	Negative
175	FFA	FA(20:4)	FALSE	303.2	Unit	303.2	Unit	3.4	2	166	0	4	Negative
176	FFA	FA(20:4) d11 (IS)	TRUE	314.2	Unit	314.2	Unit	3.4	1	166	0	4	Negative
177	FFA	FA(20:5)	FALSE	301.2	Unit	301.2	Unit	3.0	2	166	0	4	Negative
178	FFA	FA(22:4)	FALSE	331.3	Unit	331.3	Unit	3.9	2	166	0	4	Negative
179	FFA	FA(22:5)	FALSE	329.3	Unit	329.2	Unit	3.6	2	166	0	4	Negative
180	FFA	FA(22:6) d5 (IS)	TRUE	332.2	Unit	332.2	Unit	3.2	1	166	0	4	Negative
181	FFA	FA(22:6)	FALSE	327.2	Unit	327.2	Unit	3.3	2	166	0	4	Negative
182	GM3	GM3(d18:1/16:0)	FALSE	1153.7	Unit	264.3	Unit	4.9	1	166	61	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
183	GM3	GM3(d18:1/18:0)	FALSE	1181.8	Unit	264.3	Unit	5.6	1	166	61	5	Positive
184	GM3	GM3(d18:1/20:0)	FALSE	1209.8	Unit	264.3	Unit	6.5	1	166	61	5	Positive
185	GM3	GM3(d18:1/22:0)	FALSE	1237.8	Unit	264.3	Unit	7.5	1	166	61	5	Positive
186	GM3	GM3(d18:1/24:0)	FALSE	1265.8	Unit	264.3	Unit	8.5	1	166	61	5	Positive
187	GM3	GM3(d18:1/24:1)	FALSE	1263.8	Unit	264.3	Unit	7.5	1	166	61	5	Positive
188	HexCer	HexCer(d16:1/18:0)	FALSE	700.6	Unit	236.3	Unit	6.6	0.5	166	35	5	Positive
189	HexCer	HexCer(d16:1/20:0)	FALSE	728.6	Unit	236.3	Unit	7.7	0.5	166	35	5	Positive
190	HexCer	HexCer(d16:1/22:0)	FALSE	756.7	Unit	236.3	Unit	8.9	0.5	166	35	5	Positive
191	HexCer	HexCer(d16:1/24:0)	FALSE	784.7	Unit	236.3	Unit	10.0	0.5	166	35	5	Positive
192	HexCer	HexCer(d18:1/16:0)	FALSE	700.6	Unit	264.3	Unit	6.6	0.6	166	35	5	Positive
193	HexCer	HexCer(d18:1/18:0)	FALSE	728.6	Unit	264.3	Unit	7.6	0.6	166	35	5	Positive
194	HexCer	HexCer(d18:1/20:0)	FALSE	756.6	Unit	264.3	Unit	8.8	0.6	166	35	5	Positive
195	HexCer	HexCer(d18:1/22:0)	FALSE	784.7	Unit	264.3	Unit	10.0	0.6	166	35	5	Positive
196	HexCer	HexCer(d18:1/24:0)	FALSE	812.7	Unit	264.3	Unit	10.2	0.6	166	35	5	Positive
197	HexCer	HexCer(d18:1/24:1)	FALSE	810.7	Unit	264.3	Unit	10.0	0.6	166	35	5	Positive
198	HexCer	HexCer(d18:2/20:0)	FALSE	754.6	Unit	262.3	Unit	7.9	0.5	166	35	5	Positive
199	HexCer	HexCer(d18:2/22:0)	FALSE	782.7	Unit	262.3	Unit	9.0	0.5	166	35	5	Positive
200	HexCer	HexCer(d18:2/24:0)	FALSE	810.7	Unit	262.3	Unit	10.1	0.5	166	35	5	Positive
201	Hex2Cer	Hex2Cer(d16:1/16:0)	FALSE	834.6	Unit	236.3	Unit	5.2	0.5	166	49	5	Positive
202	Hex2Cer	Hex2Cer(d16:1/24:1)	FALSE	944.7	Unit	236.3	Unit	8.2	0.5	166	49	5	Positive
203	Hex2Cer	Hex2Cer(d18:1/15:0) d7 (IS)	TRUE	855.6	Unit	271.3	Unit	5.6	1	166	49	5	Positive
204	Hex2Cer	Hex2Cer(d18:1/16:0)	FALSE	862.6	Unit	264.3	Unit	6.1	0.6	166	49	5	Positive
205	Hex2Cer	Hex2Cer(d18:1/20:0)	FALSE	918.7	Unit	264.3	Unit	8.1	0.6	166	49	5	Positive
206	Hex2Cer	Hex2Cer(d18:1/22:0)	FALSE	946.7	Unit	264.3	Unit	9.3	0.6	166	49	5	Positive
207	Hex2Cer	Hex2Cer(d18:1/24:0)	FALSE	974.8	Unit	264.3	Unit	10.1	0.6	166	49	5	Positive
208	Hex2Cer	Hex2Cer(d18:1/24:1)	FALSE	972.7	Unit	264.3	Unit	9.3	0.6	166	49	5	Positive
209	Hex2Cer	Hex2Cer(d18:2/16:0)	FALSE	860.6	Unit	262.3	Unit	5.4	0.5	166	49	5	Positive
210	Hex2Cer	Hex2Cer(d18:2/24:1)	FALSE	970.7	Unit	262.3	Unit	8.4	0.5	166	49	5	Positive
211	Hex3Cer	Hex3Cer(d18:1/16:0)	FALSE	1024.7	Unit	264.3	Unit	5.7	0.5	166	61	5	Positive
212	Hex3Cer	Hex3Cer(d18:1/17:0) (IS)	TRUE	1038.7	Unit	264.3	Unit	6.2	0.5	166	61	5	Positive
213	Hex3Cer	Hex3Cer(d18:1/18:0)	FALSE	1052.7	Unit	264.3	Unit	6.6	0.5	166	61	5	Positive
214	Hex3Cer	Hex3Cer(d18:1/22:0)	FALSE	1108.8	Unit	264.3	Unit	8.8	0.5	166	61	5	Positive
215	Hex3Cer	Hex3Cer(d18:1/24:0)	FALSE	1136.8	Unit	264.3	Unit	10.0	0.7	166	61	5	Positive
216	Hex3Cer	Hex3Cer(d18:1/24:1)	FALSE	1134.8	Unit	264.3	Unit	8.8	0.5	166	61	5	Positive
217	Hex2Cer	HexCer(d18:1/15:0) d7 (IS)	TRUE	693.6	Unit	271.3	Unit	6.1	1	166	35	5	Positive
218	LPC	LPC(14:0) (a\b)	FALSE	468.3	Unit	184.1	Unit	1.9	1	166	21	5	Positive
219	LPC	LPC(15:0) (a\b)	FALSE	482.3	Unit	184.1	Unit	2.3	1	166	21	5	Positive
220	LPC	LPC(16:0) (a\b)	FALSE	496.3	Unit	184.1	Unit	2.6	1	166	21	5	Positive
221	LPC	LPC(16:1) (a\b)	FALSE	494.3	Unit	184.1	Unit	2.1	1	166	21	5	Positive
222	LPC	LPC(17:0) [sn1]	FALSE	510.4	Unit	184.1	Unit	2.9	1	166	21	5	Positive
223	LPC	LPC(17:1) (a\b\c)	FALSE	508.4	Unit	184.1	Unit	2.5	1	166	21	5	Positive
224	LPC	LPC(18:0) (a\b)	FALSE	524.4	Unit	184.1	Unit	3.3	1	166	21	5	Positive
225	LPC	LPC(18:1) (a\b)	FALSE	522.4	Unit	184.1	Unit	2.8	1	166	21	5	Positive
226	LPC	LPC(18:1) d7 (IS)	TRUE	529.4	Unit	184.1	Unit	2.8	1	166	21	5	Positive
227	LPC	LPC(18:2) (a\b)	FALSE	520.3	Unit	184.1	Unit	2.3	1	166	21	5	Positive
228	oxLipid	LPC(18:2) [+OH]	FALSE	536.3	Unit	184.1	Unit	1.0	1	166	21	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	$\Delta$ RT	碎裂电压	碰撞能量	池加速电压	极性
229	LPC	LPC(18:3) (a\b\c)	FALSE	518.3	Unit	184.1	Unit	2.0	1	166	21	5	Positive
230	LPC (104)	LPC 18:3(104)	FALSE	518.3	Unit	104.1	Unit	2.0	0.8	166	21	5	Positive
231	LPC	LPC(19:0) (a\b\c)	FALSE	538.4	Unit	184.1	Unit	3.5	1	166	21	5	Positive
232	LPC	LPC(19:1) (a\b\c)	FALSE	536.4	Unit	184.1	Unit	3.1	1	166	21	5	Positive
233	LPC	LPC(20:0) (a\b)	FALSE	552.4	Unit	184.1	Unit	3.8	1	166	21	5	Positive
234	LPC	LPC(20:1) (a\b)	FALSE	550.4	Unit	184.1	Unit	3.4	1	166	21	5	Positive
235	LPC	LPC(20:2) (a\b)	FALSE	548.4	Unit	184.1	Unit	2.9	1	166	21	5	Positive
236	LPC	LPC(20:3) (a\b)	FALSE	546.4	Unit	184.1	Unit	2.6	1	166	21	5	Positive
237	LPC (104)	LPC 20:3(104)	FALSE	546.4	Unit	104.1	Unit	2.6	1	166	21	5	Positive
238	LPC	LPC(20:4) (a\b)	FALSE	544.3	Unit	184.1	Unit	2.3	1	166	21	5	Positive
239	oxLipid	LPC(20:4) [+OH]	FALSE	560.3	Unit	184.1	Unit	1.4	1.5	166	21	5	Positive
240	LPC	LPC(20:5) (a\b)	FALSE	542.3	Unit	184.1	Unit	1.9	1	166	21	5	Positive
241	LPC	LPC(22:0) (a\b)	FALSE	580.4	Unit	184.1	Unit	4.4	1	166	21	5	Positive
242	LPC	LPC(22:1) (a\b)	FALSE	578.4	Unit	184.1	Unit	3.9	1	166	21	5	Positive
243	LPC	LPC(22:4) (a\b)	FALSE	572.4	Unit	184.1	Unit	2.8	1	166	21	5	Positive
244	LPC (104)	LPC 22:5(104)	FALSE	570.4	Unit	104.1	Unit	2.5	1.5	166	21	5	Positive
245	LPC	LPC(22:5) (a\b\c)	FALSE	570.4	Unit	184.1	Unit	2.5	1	166	21	5	Positive
246	LPC	LPC(22:6) (a\b)	FALSE	568.3	Unit	184.1	Unit	2.2	1	166	21	5	Positive
247	oxLipid	LPC(22:6) [+OH]	FALSE	584.3	Unit	184.1	Unit	1.5	1.5	166	21	5	Positive
248	LPC	LPC(24:0) (a\b)	FALSE	608.5	Unit	184.1	Unit	5.1	1	166	21	5	Positive
249	LPC	LPC(26:0) (a\b)	FALSE	636.5	Unit	184.1	Unit	5.9	1	166	21	5	Positive
250	LPCO	LPC(O-16:0)	FALSE	482.4	Unit	104.1	Unit	3.0	3	166	23	5	Positive
251	LPCO	LPC(O-18:0)	FALSE	510.4	Unit	104.1	Unit	3.6	3	166	23	5	Positive
252	LPCO	LPC(O-18:1)	FALSE	508.4	Unit	104.1	Unit	3.1	3	166	23	5	Positive
253	LPCO	LPC(O-20:0)	FALSE	538.4	Unit	104.1	Unit	4.1	3	166	23	5	Positive
254	LPCO	LPC(O-20:1)	FALSE	536.4	Unit	104.1	Unit	3.7	3	166	23	5	Positive
255	LPCO	LPC(O-22:0)	FALSE	566.5	Unit	104.1	Unit	4.8	2	166	23	5	Positive
256	LPCO	LPC(O-22:1)	FALSE	564.4	Unit	104.1	Unit	4.2	2	166	23	5	Positive
257	LPCO	LPC(O-24:0)	FALSE	594.5	Unit	104.1	Unit	5.6	1	166	23	5	Positive
258	LPCO	LPC(O-24:1)	FALSE	592.5	Unit	104.1	Unit	4.8	2	166	23	5	Positive
259	LPCO	LPC(O-24:2)	FALSE	590.5	Unit	104.1	Unit	4.3	2	166	23	5	Positive
260	LPCP	LPC(P-16:0)	FALSE	480.3	Unit	104.1	Unit	3.0	0.6	166	25	5	Positive
261	LPCP	LPC(P-17:0) (a\b)	FALSE	494.3	Unit	104.1	Unit	3.3	1	166	25	5	Positive
262	LPCP	LPC(P-18:0)	FALSE	508.3	Unit	104.1	Unit	3.6	0.6	166	25	5	Positive
263	LPCP	LPC(P-18:1)	FALSE	506.3	Unit	104.1	Unit	3.1	0.6	166	25	5	Positive
264	LPCP	LPC(P-20:0)	FALSE	536.3	Unit	104.1	Unit	4.1	0.6	166	25	5	Positive
265	LPE	LPE(16:0) (a\b)	FALSE	454.3	Unit	313.3	Unit	2.7	1	166	17	5	Positive
266	LPE	LPE(17:0) (a\b)	FALSE	468.3	Unit	327.3	Unit	3.1	1	166	17	5	Positive
267	LPE	LPE(18:0) (a\b)	FALSE	482.3	Unit	341.3	Unit	3.4	1	166	17	5	Positive
268	LPE	LPE(18:1) (a\b)	FALSE	480.3	Unit	339.3	Unit	2.9	1	166	17	5	Positive
269	LPE	LPE(18:1) d7 (IS)	TRUE	487.3	Unit	346.3	Unit	2.9	1	166	17	5	Positive
270	LPE	LPE(18:2) (a\b)	FALSE	478.3	Unit	337.3	Unit	2.4	1	166	17	5	Positive
271	LPE	LPE(20:4) (a\b)	FALSE	502.3	Unit	361.3	Unit	2.4	1	166	17	5	Positive
272	LPE	LPE(22:6) (a\b)	FALSE	526.3	Unit	385.3	Unit	2.3	1	166	17	5	Positive
273	LPEP	LPE(P-16:0)	FALSE	438.3	Unit	266.4	Unit	3.1	1	166	19	5	Positive
274	LPEP	LPE(P-18:0)	FALSE	466.3	Unit	294.4	Unit	3.7	1	166	19	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
275	LPEP	LPE(P-18:1)	FALSE	464.3	Unit	292.4	Unit	3.3	1	166	19	5	Positive
276	LPEP	LPE(P-20:0)	FALSE	494.3	Unit	322.4	Unit	4.2	1	166	19	5	Positive
277	LPI	LPI 13:0 (IS)	TRUE	548.3	Unit	271.3	Unit	1.3	1	166	21	5	Positive
278	LPI	LPI(18:0)	FALSE	618.3	Unit	341.3	Unit	2.9	1	166	21	5	Positive
279	LPI	LPI(18:1) (a\b)	FALSE	616.3	Unit	339.3	Unit	2.4	1	166	21	5	Positive
280	LPI	LPI(18:2) (a\b)	FALSE	614.3	Unit	337.3	Unit	1.9	1	166	21	5	Positive
281	LPI	LPI(20:4) (a\b)	FALSE	638.3	Unit	361.3	Unit	1.9	1	166	21	5	Positive
282	Methyl-CE	methyl-CE(18:0)	FALSE	684.6	Unit	383.3	Unit	12.2	0.4	166	10	4	Positive
283	Methyl-CE	methyl-CE(18:1)	FALSE	682.6	Unit	383.3	Unit	11.9	0.4	166	10	4	Positive
284	Methyl-CE	methyl-CE(18:2)	FALSE	680.6	Unit	383.3	Unit	11.7	0.4	166	10	4	Positive
285	Methyl-CE	methyl-CE(20:4)	FALSE	704.6	Unit	383.3	Unit	11.6	0.4	166	10	4	Positive
286	mDE	methyl-DE(18:1)	FALSE	680.6	Unit	381.4	Unit	11.7	0.6	166	10	4	Positive
287	mDE	methyl-DE(18:2)	FALSE	678.6	Unit	381.4	Unit	11.5	0.6	166	10	4	Positive
288	MG	MG(18:1) d7 (IS)	TRUE	364.2	Unit	272.2	Unit	3.7	1	166	11	4	Positive
289	PA	PA(15:0_18:1) d7 (IS)	TRUE	685.6	Unit	570.6	Unit	6.2	2	166	13	4	Positive
290	PA	PA(34:1)	FALSE	692.6	Unit	577.6	Unit	6.7	2	166	13	4	Positive
291	PA	PA(36:2)	FALSE	718.6	Unit	603.6	Unit	7.0	2	166	13	4	Positive
292	PA	PA(36:3)	FALSE	716.6	Unit	601.6	Unit	6.4	2	166	13	4	Positive
293	PA	PA(36:4)	FALSE	714.6	Unit	599.6	Unit	5.7	2	166	13	4	Positive
294	PC	PC(28:0)	FALSE	678.5	Unit	184.1	Unit	5.2	0.8	166	25	5	Positive
295	PC	PC(14:0_16:0)	FALSE	706.5	Unit	184.1	Unit	6.0	0.8	166	25	5	Positive
296	PC	PC(31:0) (a\b)	FALSE	720.6	Unit	184.1	Unit	6.4	0.8	166	25	5	Positive
297	PC	PC(31:1)	FALSE	718.5	Unit	184.1	Unit	5.8	0.6	166	25	5	Positive
298	PC	PC(16:0_16:0)	FALSE	734.6	Unit	184.1	Unit	7.0	0.6	166	25	5	Positive
299	PC	PC(32:1)	FALSE	732.6	Unit	184.1	Unit	6.2	0.6	166	25	5	Positive
300	PC	PC(32:2)	FALSE	730.5	Unit	184.1	Unit	5.6	0.6	166	25	5	Positive
301	PC	PC(33:0) (a\b)	FALSE	748.6	Unit	184.1	Unit	7.5	1.2	166	25	5	Positive
302	PC	PC(33:1)	FALSE	746.6	Unit	184.1	Unit	6.7	0.6	166	25	5	Positive
303	PC	PC(33:2)	FALSE	744.6	Unit	184.1	Unit	6.0	0.6	166	25	5	Positive
304	PC	PC(16:0_18:0)	FALSE	762.6	Unit	184.1	Unit	8.1	0.6	166	25	5	Positive
305	PC	PC(16:0_18:1)	FALSE	760.6	Unit	184.1	Unit	7.2	0.6	166	25	5	Positive
306	PC	PC(16:0_18:2)	FALSE	758.6	Unit	184.1	Unit	6.5	0.6	166	25	5	Positive
307	oxLipid	PC(34:2) [+OH]	FALSE	774.6	Unit	184.1	Unit	4.8	1	166	25	5	Positive
308	PC	PC(34:3) (a\b\c)	FALSE	756.6	Unit	184.1	Unit	5.9	1.2	166	25	5	Positive
309	PC	PC(14:0_20:4)	FALSE	754.5	Unit	184.1	Unit	5.5	0.6	166	25	5	Positive
310	PC	PC(34:5)	FALSE	752.5	Unit	184.1	Unit	5.1	0.6	166	25	5	Positive
311	PC	PC(35:1) (a\b)	FALSE	774.6	Unit	184.1	Unit	7.7	1.2	166	25	5	Positive
312	PC	PC(35:2) (a\b)	FALSE	772.6	Unit	184.1	Unit	6.9	1.2	166	25	5	Positive
313	PC	PC(35:3) (a\b\c)	FALSE	770.6	Unit	184.1	Unit	6.3	3	166	25	5	Positive
314	PC	PC(15:0_20:4)	FALSE	768.6	Unit	184.1	Unit	6.0	0.6	166	25	5	Positive
315	PC	PC(35:5)	FALSE	766.5	Unit	184.1	Unit	5.4	0.6	166	25	5	Positive
316	PC	PC(18:0_18:1)	FALSE	788.6	Unit	184.1	Unit	8.3	0.8	166	25	5	Positive
317	PC	PC(36:2) (a\b)	FALSE	786.6	Unit	184.1	Unit	7.5	1	166	25	5	Positive
318	PC	PC(36:3) (a\b\c)	FALSE	784.6	Unit	184.1	Unit	6.8	1.2	166	25	5	Positive
319	PC	PC(36:4) (a\b)	FALSE	782.6	Unit	184.1	Unit	6.3	1.6	166	25	5	Positive
320	oxLipid	PC(36:4) [+OH]	FALSE	798.6	Unit	184.1	Unit	5.4	2	166	25	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
321	PC	PC(36:5) (a\b)	FALSE	780.6	Unit	184.1	Unit	5.8	1	166	25	5	Positive
322	PC	PC(36:6) (a\b)	FALSE	778.5	Unit	184.1	Unit	5.3	0.6	166	25	5	Positive
323	PC	PC(37:4) (a\b)	FALSE	796.6	Unit	184.1	Unit	6.9	1.2	166	25	5	Positive
324	PC	PC(15:0_22:6)	FALSE	792.6	Unit	184.1	Unit	5.8	0.6	166	25	5	Positive
325	PC	PC(38:2)	FALSE	814.6	Unit	184.1	Unit	8.6	1.4	166	25	5	Positive
326	PC	PC(18:0_20:3)	FALSE	812.6	Unit	184.1	Unit	7.9	1.4	166	25	5	Positive
327	PC	PC(38:4) (a\b\c)	FALSE	810.6	Unit	184.1	Unit	7.3	1.6	166	25	5	Positive
328	PC	PC(38:5) (a\b)	FALSE	808.6	Unit	184.1	Unit	6.7	1	166	25	5	Positive
329	PC	PC(38:6) (a\b)	FALSE	806.6	Unit	184.1	Unit	6.1	1.2	166	25	5	Positive
330	oxLipid	PC(38:6) [+OH]	FALSE	822.6	Unit	184.1	Unit	5.3	2	166	25	5	Positive
331	PC	PC(38:7) (a\b\c)	FALSE	804.6	Unit	184.1	Unit	5.6	1	166	25	5	Positive
332	PC	PC(39:5) (a\b)	FALSE	822.6	Unit	184.1	Unit	6.8	1.2	166	25	5	Positive
333	PC	PC(39:6) (a\b)	FALSE	820.6	Unit	184.1	Unit	6.7	1	166	25	5	Positive
334	PC	PC(40:4) (a\b)	FALSE	838.6	Unit	184.1	Unit	8.4	1.2	166	25	5	Positive
335	PC	PC(40:5) (a\b)	FALSE	836.6	Unit	184.1	Unit	7.8	1.2	166	25	5	Positive
336	PC	PC(18:0_22:6)	FALSE	834.6	Unit	184.1	Unit	7.3	0.6	166	25	5	Positive
337	PC	PC(40:7) (a\b\c)	FALSE	832.6	Unit	184.1	Unit	6.4	1	166	25	5	Positive
338	PC	PC(40:8)	FALSE	830.6	Unit	184.1	Unit	5.8	0.8	166	25	5	Positive
339	PC	PC(15:0_18:1) d7 (IS)	TRUE	753.6	Unit	184.1	Unit	6.7	0.7	166	25	5	Positive
340	PC (O)	PC(O-16:0/16:0)	FALSE	720.6	Unit	184.1	Unit	7.7	0.8	166	25	5	Positive
341	PC (O)	PC(O-32:1)	FALSE	718.5	Unit	184.1	Unit	7.2	0.8	166	25	5	Positive
342	PC (O)	PC(O-32:2)	FALSE	716.6	Unit	184.1	Unit	6.3	0.8	166	25	5	Positive
343	PC (O)	PC(O-34:1)	FALSE	746.6	Unit	184.1	Unit	7.9	0.8	166	25	5	Positive
344	PC (O)	PC(O-34:2)	FALSE	744.6	Unit	184.1	Unit	7.2	1	166	25	5	Positive
345	PC (O)	PC(O-34:4)	FALSE	740.6	Unit	184.1	Unit	6.1	0.8	166	25	5	Positive
346	PC (O)	PC(O-35:4)	FALSE	754.5	Unit	184.1	Unit	6.7	0.8	166	25	5	Positive
347	PC (O)	PC(O-36:0)	FALSE	776.6	Unit	184.1	Unit	10.0	0.8	166	25	5	Positive
348	PC (O)	PC(O-18:0/18:1)	FALSE	774.6	Unit	184.1	Unit	9.1	0.8	166	25	5	Positive
349	PC (O)	PC(O-36:2) (a\b)	FALSE	772.6	Unit	184.1	Unit	8.2	1.2	166	25	5	Positive
350	PC (O)	PC(O-36:3) (a\b)	FALSE	770.6	Unit	184.1	Unit	7.4	1.2	166	25	5	Positive
351	PC (O)	PC(O-16:0/20:4)	FALSE	768.6	Unit	184.1	Unit	7.1	0.8	166	25	5	Positive
352	PC (O)	PC(O-36:5)	FALSE	766.5	Unit	184.1	Unit	6.4	0.8	166	25	5	Positive
353	PC (O)	PC(O-18:0/20:4)	FALSE	796.6	Unit	184.1	Unit	8.2	0.8	166	25	5	Positive
354	PC (O)	PC(O-38:5)	FALSE	794.6	Unit	184.1	Unit	7.2	0.8	166	25	5	Positive
355	PC (O)	PC(O-16:0/22:6)	FALSE	792.6	Unit	184.1	Unit	6.9	0.8	166	25	5	Positive
356	PC (O)	PC(O-40:5)	FALSE	822.6	Unit	184.1	Unit	8.3	0.8	166	25	5	Positive
357	PC (O)	PC(O-18:0/22:6)	FALSE	820.6	Unit	184.1	Unit	8.0	1	166	25	5	Positive
358	PC (O)	PC(O-40:7)	FALSE	818.6	Unit	184.1	Unit	7.0	1	166	25	5	Positive
359	PC (P)	PC(P-18:0/18:1) d9 (IS)	TRUE	781.6	Unit	184.1	Unit	8.9	1	166	25	5	Positive
360	PC (P)	PC(P-16:0/14:0)	FALSE	690.4	Unit	184.1	Unit	6.6	0.8	166	25	5	Positive
361	PC (P)	PC(P-16:0/16:0)	FALSE	718.5	Unit	184.1	Unit	7.6	0.8	166	25	5	Positive
362	PC (P)	PC(P-16:0/16:1)	FALSE	716.6	Unit	184.1	Unit	6.8	0.8	166	25	5	Positive
363	PC (P)	PC(P-16:0/18:0)	FALSE	746.6	Unit	184.1	Unit	8.8	0.8	166	25	5	Positive
364	PC (P)	PC(P-16:0/18:1)	FALSE	744.6	Unit	184.1	Unit	7.8	0.8	166	25	5	Positive
365	PC (P)	PC(P-16:0/18:2)	FALSE	742.5	Unit	184.1	Unit	7.0	0.8	166	25	5	Positive
366	PC (P)	PC(P-16:0/18:3)	FALSE	740.6	Unit	184.1	Unit	6.4	0.8	166	25	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
367	PC (P)	PC(P-35:2) (a\b)	FALSE	756.6	Unit	184.1	Unit	7.5	1.5	166	25	5	Positive
368	PC (P)	PC(P-15:0/20:4) (a\b)	FALSE	752.6	Unit	184.1	Unit	6.3	1	166	25	5	Positive
369	PC (P)	PC(P-36:2) (a\b)	FALSE	770.6	Unit	184.1	Unit	8.1	1	166	25	5	Positive
370	PC (P)	PC(P-36:3)	FALSE	768.5	Unit	184.1	Unit	7.3	1	166	25	5	Positive
371	PC (P)	PC(P-16:0/20:4)	FALSE	766.5	Unit	184.1	Unit	6.9	0.8	166	25	5	Positive
372	PC (P)	PC(P-16:0/20:5)	FALSE	764.6	Unit	184.1	Unit	6.3	0.8	166	25	5	Positive
373	PC (P)	PC(P-37:4) (a\b)	FALSE	780.5	Unit	184.1	Unit	7.4	1	166	25	5	Positive
374	PC (P)	PC(P-18:0/20:4)	FALSE	794.6	Unit	184.1	Unit	8.0	0.8	166	25	5	Positive
375	PC (P)	PC(P-38:5) (a\b)	FALSE	792.6	Unit	184.1	Unit	7.2	1.2	166	25	5	Positive
376	PC (P)	PC(P-16:0/22:6)	FALSE	790.6	Unit	184.1	Unit	6.7	0.8	166	25	5	Positive
377	PC (P)	PC(P-20:0/20:4)	FALSE	822.6	Unit	184.1	Unit	9.2	0.8	166	25	5	Positive
378	PC (P)	PC(P-18:0/22:5)	FALSE	820.6	Unit	184.1	Unit	8.1	1	166	25	5	Positive
379	PC (P)	PC(P-18:0/22:6)	FALSE	818.6	Unit	184.1	Unit	7.8	0.8	166	25	5	Positive
380	PC (P)	PC(P-18:1/22:6)	FALSE	816.6	Unit	184.1	Unit	6.8	0.8	166	25	5	Positive
381	PE	PE(16:0_16:0)	FALSE	692.5	Unit	551.5	Unit	7.4	0.6	166	19	5	Positive
382	PE	PE(16:0_16:1)	FALSE	690.5	Unit	549.5	Unit	6.6	0.6	166	19	5	Positive
383	PE	PE(16:0_18:1)	FALSE	718.5	Unit	577.5	Unit	7.6	0.6	166	19	5	Positive
384	PE	PE(16:0_18:2)	FALSE	716.5	Unit	575.5	Unit	6.8	0.6	166	19	5	Positive
385	PE	PE(34:3) (a\b\c)	FALSE	714.5	Unit	573.5	Unit	6.2	1	166	19	5	Positive
386	PE	PE(35:1) (a\b)	FALSE	732.6	Unit	591.5	Unit	8.0	1	166	19	5	Positive
387	PE	PE(35:2) (a\b)	FALSE	730.5	Unit	589.5	Unit	7.3	1	166	19	5	Positive
388	PE	PE(18:0_18:1)	FALSE	746.6	Unit	605.6	Unit	8.8	0.6	166	19	5	Positive
389	PE	PE(36:2) (a\b)	FALSE	744.6	Unit	603.5	Unit	7.9	1	166	19	5	Positive
390	PE	PE(36:3) (a\b)	FALSE	742.5	Unit	601.5	Unit	7.1	0.6	166	19	5	Positive
391	PE	PE(16:0_20:4)	FALSE	740.5	Unit	599.5	Unit	6.8	0.6	166	19	5	Positive
392	PE	PE(36:5) (a\b)	FALSE	738.5	Unit	597.5	Unit	6.1	1	166	19	5	Positive
393	PE	PE(37:4) (a\b)	FALSE	754.6	Unit	613.5	Unit	7.2	1	166	19	5	Positive
394	PE	PE(18:0_20:3) (a\b)	FALSE	770.6	Unit	629.6	Unit	8.5	1.2	166	19	5	Positive
395	PE	PE(18:0_20:4)	FALSE	768.6	Unit	627.5	Unit	7.9	0.6	166	19	5	Positive
396	PE	PE(38:5) (a\b)	FALSE	766.5	Unit	625.5	Unit	7.1	1	166	19	5	Positive
397	PE	PE(16:0_22:6)	FALSE	764.5	Unit	623.5	Unit	6.6	0.6	166	19	5	Positive
398	PE	PE(39:6) (a\b)	FALSE	778.5	Unit	637.5	Unit	7.0	1	166	19	5	Positive
399	PE	PE(40:4) (a\b)	FALSE	796.6	Unit	655.6	Unit	8.9	1	166	19	5	Positive
400	PE	PE(18:0_22:5) (a\b)	FALSE	794.6	Unit	653.6	Unit	8.2	1.4	166	19	5	Positive
401	PE	PE(18:0_22:6)	FALSE	792.6	Unit	651.5	Unit	7.6	0.6	166	19	5	Positive
402	PE	PE(18:1_22:6) (a\b)	FALSE	790.5	Unit	649.5	Unit	6.8	0.6	166	19	5	Positive
403	PE	PE(15:0_18:1) d7 (IS)	TRUE	711.6	Unit	570.5	Unit	7.0	1	166	19	5	Positive
404	PE (O)	PE(O-34:1)	FALSE	704.6	Unit	563.5	Unit	8.3	0.8	166	19	5	Positive
405	PE (O)	PE(O-16:0/18:2)	FALSE	702.5	Unit	561.5	Unit	7.6	0.8	166	19	5	Positive
406	PE (O)	PE(O-36:3) (a\b)	FALSE	728.6	Unit	587.5	Unit	7.8	1	166	19	5	Positive
407	PE (O)	PE(O-16:0/20:4)	FALSE	726.5	Unit	585.5	Unit	7.5	0.8	166	19	5	Positive
408	PE (O)	PE(O-36:5)	FALSE	724.5	Unit	583.5	Unit	6.8	0.8	166	19	5	Positive
409	PE (O)	PE(O-38:4) (a\b)	FALSE	754.6	Unit	613.6	Unit	8.5	1.2	166	19	5	Positive
410	PE (O)	PE(O-38:5) (a\b)	FALSE	752.6	Unit	611.5	Unit	7.8	1.2	166	19	5	Positive
411	PE (O)	PE(O-16:0/22:6)	FALSE	750.6	Unit	609.5	Unit	7.3	1.2	166	19	5	Positive
412	PE (O)	PE(O-18:0/22:5)	FALSE	780.6	Unit	639.6	Unit	8.7	0.8	166	19	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
413	PE (O)	PE(O-18:0/22:6)	FALSE	778.5	Unit	637.5	Unit	8.2	0.8	166	19	5	Positive
414	PE (O)	PE(O-18:1/22:6)	FALSE	776.6	Unit	635.5	Unit	7.4	0.8	166	19	5	Positive
415	PE (P)	PE(P-15:0/22:6)	FALSE	734.5	Unit	385.3	Unit	6.5	1	166	19	5	Positive
416	PE (P)	PE(P-16:0/18:1)	FALSE	702.5	Unit	339.3	Unit	8.3	0.6	166	19	5	Positive
417	PE (P)	PE(P-16:0/18:2)	FALSE	700.5	Unit	337.3	Unit	7.4	0.6	166	19	5	Positive
418	PE (P)	PE(P-16:0/20:3) (a\b)	FALSE	726.5	Unit	363.3	Unit	7.8	0.6	166	19	5	Positive
419	PE (P)	PE(P-16:0/20:4)	FALSE	724.5	Unit	361.3	Unit	7.3	0.6	166	19	5	Positive
420	PE (P)	PE(P-16:0/20:5)	FALSE	722.5	Unit	359.3	Unit	6.7	0.6	166	19	5	Positive
421	PE (P)	PE(P-16:0/22:4)	FALSE	752.6	Unit	389.3	Unit	8.2	0.6	166	19	5	Positive
422	PE (P)	PE(P-16:0/22:5) (a\b)	FALSE	750.5	Unit	387.3	Unit	7.6	1.4	166	19	5	Positive
423	PE (P)	PE(P-16:0/22:6)	FALSE	748.5	Unit	385.3	Unit	7.1	0.6	166	19	5	Positive
424	PE (P)	PE(P-17:0/20:4) (a\b)	FALSE	738.6	Unit	361.3	Unit	7.8	1	166	19	5	Positive
425	PE (P)	PE(P-17:0/22:6) (a\b)	FALSE	762.6	Unit	385.3	Unit	7.6	1	166	19	5	Positive
426	PE (P)	PE(P-18:0/18:1)	FALSE	730.6	Unit	339.3	Unit	9.5	0.6	166	19	5	Positive
427	PE (P)	PE(P-18:0/18:1) d9 (IS)	TRUE	739.5	Unit	348.3	Unit	9.4	1	166	19	5	Positive
428	PE (P)	PE(P-18:0/18:2)	FALSE	728.6	Unit	337.3	Unit	8.6	0.6	166	19	5	Positive
429	PE (P)	PE(P-18:0/18:3)	FALSE	726.5	Unit	335.3	Unit	7.9	1	166	19	5	Positive
430	PE (P)	PE(P-18:0/20:3) (a\b)	FALSE	754.5	Unit	363.3	Unit	9.0	0.6	166	19	5	Positive
431	PE (P)	PE(P-18:0/20:4)	FALSE	752.6	Unit	361.3	Unit	8.5	0.6	166	19	5	Positive
432	PE (P)	PE(P-18:0/20:5)	FALSE	750.5	Unit	359.3	Unit	7.8	0.6	166	19	5	Positive
433	PE (P)	PE(P-18:0/22:4)	FALSE	780.6	Unit	389.3	Unit	9.4	0.6	166	19	5	Positive
434	PE (P)	PE(P-18:0/22:5) (a\b)	FALSE	778.5	Unit	387.3	Unit	8.8	1.4	166	19	5	Positive
435	PE (P)	PE(P-18:0/22:6)	FALSE	776.6	Unit	385.3	Unit	8.2	0.6	166	19	5	Positive
436	PE (P)	PE(P-18:1/18:1) (a\b)	FALSE	728.6	Unit	339.3	Unit	8.5	0.6	166	19	5	Positive
437	PE (P)	PE(P-18:1/18:2) (a\b)	FALSE	726.5	Unit	337.3	Unit	7.7	0.6	166	19	5	Positive
438	PE (P)	PE(P-18:1/20:3) (a\b)	FALSE	752.5	Unit	363.3	Unit	8.0	0.6	166	19	5	Positive
439	PE (P)	PE(P-18:1/20:4) (a\b)	FALSE	750.5	Unit	361.3	Unit	7.6	0.6	166	19	5	Positive
440	PE (P)	PE(P-18:1/20:5) (a\b)	FALSE	748.5	Unit	359.3	Unit	6.9	0.6	166	19	5	Positive
441	PE (P)	PE(P-18:1/22:4)	FALSE	778.5	Unit	389.3	Unit	8.4	0.6	166	19	5	Positive
442	PE (P)	PE(P-18:1/22:5) (a\b)	FALSE	776.6	Unit	387.3	Unit	7.8	1.4	166	19	5	Positive
443	PE (P)	PE(P-18:1/22:6) (a\b)	FALSE	774.5	Unit	385.3	Unit	7.3	1	166	19	5	Positive
444	PE (P)	PE(P-19:0/20:4) (a\b)	FALSE	766.6	Unit	361.3	Unit	9.0	1	166	19	5	Positive
445	PE (P)	PE(P-20:0/18:1)	FALSE	758.6	Unit	339.3	Unit	10.1	0.6	166	19	5	Positive
446	PE (P)	PE(P-20:0/18:2)	FALSE	756.6	Unit	337.3	Unit	9.9	0.6	166	19	5	Positive
447	PE (P)	PE(P-20:0/20:4)	FALSE	780.6	Unit	361.3	Unit	9.7	0.6	166	19	5	Positive
448	PE (P)	PE(P-20:0/22:6)	FALSE	804.6	Unit	385.3	Unit	9.5	0.6	166	19	5	Positive
449	PE (P)	PE(P-20:1/20:4)	FALSE	778.5	Unit	361.3	Unit	8.6	0.6	166	19	5	Positive
450	PE (P)	PE(P-20:1/22:6)	FALSE	802.6	Unit	385.3	Unit	8.3	0.6	166	19	5	Positive
451	PG	PG(34:1)	FALSE	766.6	Unit	577.5	Unit	6.0	0.6	166	17	5	Positive
452	PG	PG(36:1)	FALSE	794.6	Unit	605.6	Unit	7.1	0.6	166	17	5	Positive
453	PG	PG(36:2)	FALSE	792.6	Unit	603.5	Unit	6.5	0.6	166	17	5	Positive
454	PG	PG(15:0_18:1) d7 (IS)	TRUE	759.6	Unit	570.6	Unit	5.9	1	166	17	5	Positive
455	PI	PI(15:0_18:1) d7 (IS)	FALSE	847.6	Unit	570.6	Unit	5.6	1.5	166	17	5	Positive
456	PI	PI(16:0/16:0)	FALSE	828.6	Unit	551.6	Unit	5.9	0.8	166	17	5	Positive
457	PI	PI(16:0_16:1)	FALSE	826.5	Unit	549.5	Unit	5.3	0.8	166	17	5	Positive
458	PI	PI(34:0)	FALSE	856.6	Unit	579.6	Unit	6.8	0.8	166	17	5	Positive



化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
459	PI	PI(34:1)	FALSE	854.6	Unit	577.6	Unit	6.0	0.8	166	17	5	Positive
460	PI	PI(15-MHDA_18:1)/PI(17:0_18:1)	FALSE	868.6	Unit	591.6	Unit	6.5	0.8	166	17	5	Positive
461	PI	PI(15-MHDA_18:2)/PI(17:0_18:2)	FALSE	866.6	Unit	589.6	Unit	5.9	0.8	166	17	5	Positive
462	PI	PI(18:0_18:1)	FALSE	882.6	Unit	605.6	Unit	7.0	0.8	166	17	5	Positive
463	PI	PI(36:2)	FALSE	880.6	Unit	603.6	Unit	6.3	3	166	17	5	Positive
464	PI	PI(36:3) (a\b)	FALSE	878.6	Unit	601.6	Unit	5.7	1	166	17	5	Positive
465	PI	PI(16:0_20:4)	FALSE	876.6	Unit	599.6	Unit	5.4	0.8	166	17	5	Positive
466	PI	PI(37:4)	FALSE	890.6	Unit	613.6	Unit	5.8	1	166	17	5	Positive
467	PI	PI(37:6)	FALSE	886.6	Unit	609.6	Unit	5.7	1	166	17	5	Positive
468	PI	PI(18:0_20:2)	FALSE	908.6	Unit	631.6	Unit	7.2	0.8	166	17	5	Positive
469	PI	PI(18:0_20:3) (a\b)	FALSE	906.6	Unit	629.6	Unit	6.8	1	166	17	5	Positive
470	PI	PI(18:0_20:4)	FALSE	904.6	Unit	627.6	Unit	6.3	0.8	166	17	5	Positive
471	PI	PI(38:5) (a\b)	FALSE	902.6	Unit	625.6	Unit	5.6	1	166	17	5	Positive
472	PI	PI(38:6)	FALSE	900.6	Unit	623.6	Unit	5.3	0.8	166	17	5	Positive
473	PI	PI(40:4) (a\b)	FALSE	932.6	Unit	655.6	Unit	7.1	0.8	166	17	5	Positive
474	PI	PI(18:0_22:5) (a\b)	FALSE	930.6	Unit	653.6	Unit	6.5	1	166	17	5	Positive
475	PI	PI(18:0_22:6)	FALSE	928.6	Unit	651.6	Unit	6.1	0.8	166	17	5	Positive
476	PIP	PIP1(38:4)	FALSE	984.7	Unit	627.7	Unit	5.3	2	166	20	4	Positive
477	PS	PS(36:1)	FALSE	790.6	Unit	605.6	Unit	7.1	1	166	23	5	Positive
478	PS	PS(36:2)	FALSE	788.5	Unit	603.5	Unit	6.4	1	166	23	5	Positive
479	PS	PS(38:4)	FALSE	812.5	Unit	627.5	Unit	6.3	1	166	23	5	Positive
480	PS	PS(40:5)	FALSE	838.6	Unit	653.6	Unit	6.4	1	166	23	5	Positive
481	PS	PS(15:0_18:1) d7 (IS)	TRUE	755.5	Unit	570.5	Unit	5.7	1	166	23	4	Positive
482	S1P	S1P(d16:1)	FALSE	352.2	Unit	236.3	Unit	1.5	3	166	12	4	Positive
483	S1P	S1P(d18:0)	FALSE	382.2	Unit	284.3	Unit	2.3	3	166	12	4	Positive
484	S1P	S1P(d18:1)	FALSE	380.2	Unit	264.3	Unit	2.1	3	166	12	4	Positive
485	S1P	S1P(d18:1) d7 (IS)	TRUE	387.2	Unit	271.3	Unit	2.1	1	166	12	4	Positive
486	S1P	S1P(d18:2)	FALSE	378.2	Unit	262.3	Unit	1.6	3	166	12	4	Positive
487	SM	SM(d17:1/14:0)	FALSE	661.5	Unit	184.1	Unit	4.8	0.6	166	29	5	Positive
488	SM	SM(d18:0/14:0)	FALSE	677.6	Unit	184.1	Unit	5.5	0.6	166	29	5	Positive
489	SM	SM(d18:1/14:0)/SM(d16:1/16:0)	FALSE	675.5	Unit	184.1	Unit	5.2	0.6	166	29	5	Positive
490	SM	SM(d18:2/14:0)	FALSE	673.5	Unit	184.1	Unit	4.7	0.6	166	29	5	Positive
491	SM	SM(d17:1/16:0)	FALSE	689.6	Unit	184.1	Unit	5.6	0.6	166	29	5	Positive
492	SM	SM(d18:0/16:0)	FALSE	705.6	Unit	184.1	Unit	6.4	0.6	166	29	5	Positive
493	SM	SM(d18:1/16:0)	FALSE	703.6	Unit	184.1	Unit	6.0	0.6	166	29	5	Positive
494	SM	SM(d18:2/16:0)	FALSE	701.6	Unit	184.1	Unit	5.4	0.6	166	29	5	Positive
495	SM	SM(34:3)	FALSE	699.5	Unit	184.1	Unit	4.9	0.6	166	29	5	Positive
496	SM	SM(35:1) (a\b)	FALSE	717.6	Unit	184.1	Unit	6.4	0.6	166	29	5	Positive
497	SM	SM(35:2) (a\b)	FALSE	715.6	Unit	184.1	Unit	5.9	1	166	29	5	Positive
498	SM	SM(d18:1/18:0)/SM(d16:1/20:0)	FALSE	731.6	Unit	184.1	Unit	7.0	0.6	166	29	5	Positive
499	SM	SM(d18:2/18:0)	FALSE	729.6	Unit	184.1	Unit	6.2	0.6	166	29	5	Positive
500	SM	SM(d18:2/18:1)	FALSE	727.6	Unit	184.1	Unit	5.6	0.6	166	29	5	Positive
501	SM	SM(37:1)	FALSE	745.6	Unit	184.1	Unit	7.6	1	166	29	5	Positive
502	SM	SM(37:2)	FALSE	743.5	Unit	184.1	Unit	6.7	0.6	166	29	5	Positive
503	SM	SM(d18:1/20:0)/SM(d16:1/22:0)	FALSE	759.6	Unit	184.1	Unit	8.2	0.6	166	29	5	Positive
504	SM	SM(d18:2/20:0)	FALSE	757.6	Unit	184.1	Unit	7.3	0.6	166	29	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
505	SM	SM(38:3) (a\b)	FALSE	755.6	Unit	184.1	Unit	6.5	1	166	29	5	Positive
506	SM	SM(d16:1/23:0)/SM(d17:1/22:0)	FALSE	773.7	Unit	184.1	Unit	8.8	0.6	166	29	5	Positive
507	SM	SM(d18:0/22:0)	FALSE	789.7	Unit	184.1	Unit	9.8	0.6	166	29	5	Positive
508	SM	SM(d18:1/22:0)/SM(d16:1/24:0)	FALSE	787.7	Unit	184.1	Unit	9.4	0.6	166	29	5	Positive
509	SM	SM(40:2) (a\b)	FALSE	785.7	Unit	184.1	Unit	8.3	1.4	166	29	5	Positive
510	SM	SM(40:3) (a\b)	FALSE	783.6	Unit	184.1	Unit	7.5	1.2	166	29	5	Positive
511	SM	SM(41:0)	FALSE	803.7	Unit	184.1	Unit	10.1	0.8	166	29	5	Positive
512	SM	SM(41:1) (a\b)	FALSE	801.7	Unit	184.1	Unit	9.9	1	166	29	5	Positive
513	SM	SM(41:2) (a\b)	FALSE	799.7	Unit	184.1	Unit	8.9	1.4	166	29	5	Positive
514	SM	SM(d18:1/24:0)	FALSE	815.7	Unit	184.1	Unit	10.1	0.6	166	29	5	Positive
515	SM	SM(42:2) (a\b)	FALSE	813.7	Unit	184.1	Unit	9.5	1.4	166	29	5	Positive
516	SM	SM(43:1)	FALSE	829.7	Unit	184.1	Unit	10.2	0.8	166	29	5	Positive
517	SM	SM(43:2) (a\b\c)	FALSE	827.7	Unit	184.1	Unit	9.9	1.4	166	29	5	Positive
518	SM	SM(44:1)	FALSE	843.6	Unit	184.1	Unit	10.3	1	166	29	5	Positive
519	SM	SM(44:2)	FALSE	841.6	Unit	184.1	Unit	10.1	1	166	29	5	Positive
520	SM	SM(44:3)	FALSE	839.6	Unit	184.1	Unit	9.6	2	166	29	5	Positive
521	SM	SM(d18:1/15:0) d9 (IS)	TRUE	698.6	Unit	193.1	Unit	5.6	1	166	29	5	Positive
522	SPN	Sph(d17:1) (IS)	FALSE	286.3	Unit	268.3	Unit	1.8	3	166	10	4	Positive
523	SPN	Sph(d18:1)	FALSE	300.3	Unit	282.3	Unit	2.4	3	166	10	4	Positive
524	SPN	Sph(d18:2)	FALSE	298.3	Unit	280.3	Unit	1.8	3	166	10	4	Positive
525	Sulfatides	SHexCer(d18:1/12:0) (IS)	TRUE	724.8	Unit	264.3	Unit	4.1	0.6	166	56	5	Positive
526	Sulfatides	SHexCer(d18:1/16:0(OH))	FALSE	796.8	Unit	264.3	Unit	5.0	0.6	166	56	5	Positive
527	Sulfatides	SHexCer(d18:1/16:0)	FALSE	780.8	Unit	264.3	Unit	5.2	0.6	166	56	5	Positive
528	Sulfatides	SHexCer(d18:1/24:0(OH))	FALSE	908.8	Unit	264.3	Unit	8.6	0.6	166	56	5	Positive
529	Sulfatides	SHexCer(d18:1/24:1(OH))	FALSE	906.8	Unit	264.3	Unit	7.5	0.8	166	56	5	Positive
530	Sulfatides	SHexCer(d18:1/24:1)	FALSE	890.8	Unit	264.3	Unit	7.7	0.6	166	56	5	Positive
531	TG	TG(48:0) [NL-16:0]	FALSE	824.8	Unit	551.5	Unit	11.4	0.7	166	25	5	Positive
532	TG	TG(48:0) [NL-18:0]	FALSE	824.8	Unit	523.5	Unit	11.4	0.7	166	25	5	Positive
533	TG	TG(48:1) [NL-16:1]	FALSE	822.8	Unit	551.5	Unit	11.3	0.7	166	25	5	Positive
534	TG	TG(48:1) [NL-18:1]	FALSE	822.8	Unit	523.5	Unit	11.3	0.7	166	25	5	Positive
535	TG	TG(48:1) [NL-18:1] d7 (IS)	TRUE	829.8	Unit	523.5	Unit	11.3	1	166	25	5	Positive
536	TG	TG(48:2) [NL-14:0]	FALSE	820.8	Unit	575.5	Unit	11.1	0.7	166	25	5	Positive
537	TG	TG(48:2) [NL-14:1]	FALSE	820.8	Unit	577.5	Unit	11.1	0.7	166	25	5	Positive
538	TG	TG(48:2) [NL-16:1]	FALSE	820.8	Unit	549.5	Unit	11.1	0.7	166	25	5	Positive
539	TG	TG(48:2) [NL-18:2]	FALSE	820.8	Unit	523.5	Unit	11.1	0.7	166	25	5	Positive
540	TG	TG(48:3) [NL-14:0]	FALSE	818.8	Unit	573.5	Unit	11.0	0.7	166	25	5	Positive
541	TG	TG(48:3) [NL-16:1]	FALSE	818.8	Unit	547.5	Unit	11.0	0.7	166	25	5	Positive
542	TG	TG(48:3) [NL-18:3]	FALSE	818.8	Unit	523.5	Unit	11.0	0.7	166	25	5	Positive
543	TG	TG(49:1) [NL-16:1]	FALSE	836.8	Unit	565.5	Unit	11.4	0.7	166	25	5	Positive
544	TG	TG(49:1) [NL-17:1]	FALSE	836.8	Unit	551.5	Unit	11.4	0.7	166	25	5	Positive
545	TG	TG(50:0) [NL-18:0]	FALSE	852.8	Unit	551.5	Unit	11.6	0.7	166	25	5	Positive
546	TG	TG(50:1) [NL-14:0]	FALSE	850.8	Unit	605.5	Unit	11.5	0.7	166	25	5	Positive
547	TG	TG(50:1) [NL-16:0]	FALSE	850.8	Unit	577.5	Unit	11.5	0.7	166	25	5	Positive
548	TG	TG(50:1) [NL-18:1]	FALSE	850.8	Unit	551.5	Unit	11.5	0.7	166	25	5	Positive
549	TG	TG(50:2) [NL-14:0]	FALSE	848.8	Unit	603.5	Unit	11.3	0.7	166	25	5	Positive
550	TG	TG(50:2) [NL-16:1]	FALSE	848.8	Unit	577.5	Unit	11.3	0.7	166	25	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	$\Delta RT$	碎裂电压	碰撞能量	池加速电压	极性
551	TG	TG(50:2) [NL-18:1]	FALSE	848.8	Unit	549.5	Unit	11.3	0.7	166	25	5	Positive
552	TG	TG(50:2) [NL-18:2]	FALSE	848.8	Unit	551.5	Unit	11.3	0.7	166	25	5	Positive
553	TG	TG(50:3) [NL-14:0]	FALSE	846.8	Unit	601.5	Unit	11.1	0.7	166	25	5	Positive
554	TG	TG(50:3) [NL-14:1]	FALSE	846.8	Unit	603.5	Unit	11.1	0.7	166	25	5	Positive
555	TG	TG(50:3) [NL-16:1]	FALSE	846.8	Unit	575.5	Unit	11.1	0.7	166	25	5	Positive
556	TG	TG(50:3) [NL-18:2]	FALSE	846.8	Unit	549.5	Unit	11.1	0.7	166	25	5	Positive
557	TG	TG(50:3) [NL-18:3]	FALSE	846.8	Unit	551.5	Unit	11.2	0.7	166	25	5	Positive
558	TG	TG(50:4) [NL-14:0]	FALSE	844.8	Unit	599.5	Unit	11.0	0.7	166	25	5	Positive
559	TG	TG(50:4) [NL-18:3]	FALSE	844.8	Unit	549.5	Unit	11.0	0.7	166	25	5	Positive
560	TG	TG(50:4) [NL-20:4]	FALSE	844.8	Unit	523.5	Unit	11.1	0.7	166	25	5	Positive
561	TG	TG(51:0) [NL-16:0]	FALSE	866.7	Unit	593.4	Unit	11.7	0.7	166	25	5	Positive
562	TG	TG(51:1) [NL-17:0]	FALSE	864.8	Unit	577.5	Unit	11.5	0.7	166	25	5	Positive
563	TG	TG(51:2) [NL-15:0]	FALSE	862.8	Unit	603.5	Unit	11.4	0.7	166	25	5	Positive
564	TG	TG(51:2) [NL-17:0]	FALSE	862.8	Unit	575.5	Unit	11.4	0.7	166	25	5	Positive
565	TG	TG(51:2) [NL-17:1]	FALSE	862.8	Unit	577.5	Unit	11.4	0.7	166	25	5	Positive
566	TG	TG(52:1) [NL-18:0]	FALSE	878.8	Unit	577.5	Unit	11.7	0.7	166	25	5	Positive
567	TG	TG(52:1) [NL-18:1]	FALSE	878.8	Unit	579.5	Unit	11.7	0.7	166	25	5	Positive
568	TG	TG(52:2) [NL-16:0]	FALSE	876.8	Unit	603.5	Unit	11.5	0.7	166	25	5	Positive
569	TG	TG(52:2) [NL-18:2]	FALSE	876.8	Unit	579.5	Unit	11.5	0.7	166	25	5	Positive
570	TG	TG(52:3) [NL-16:1]	FALSE	874.8	Unit	603.5	Unit	11.3	0.7	166	25	5	Positive
571	TG	TG(52:3) [NL-18:2]	FALSE	874.8	Unit	577.5	Unit	11.3	0.7	166	25	5	Positive
572	TG	TG(52:4) [NL-16:1]	FALSE	872.8	Unit	601.5	Unit	11.2	0.7	166	25	5	Positive
573	TG	TG(52:4) [NL-18:2]	FALSE	872.8	Unit	575.5	Unit	11.2	0.7	166	25	5	Positive
574	TG	TG(52:4) [NL-18:3]	FALSE	872.8	Unit	577.5	Unit	11.2	0.7	166	25	5	Positive
575	TG	TG(52:5) [NL-18:3]	FALSE	870.8	Unit	575.5	Unit	11.1	0.7	166	25	5	Positive
576	TG	TG(52:5) [NL-20:4]	FALSE	870.8	Unit	549.5	Unit	11.1	0.7	166	25	5	Positive
577	TG	TG(52:5) [NL-20:5]	FALSE	870.8	Unit	551.5	Unit	11.1	0.7	166	25	5	Positive
578	TG	TG(53:2) [NL-17:1]	FALSE	890.8	Unit	605.5	Unit	11.6	0.7	166	25	5	Positive
579	TG	TG(53:2) [NL-18:1]	FALSE	890.8	Unit	591.5	Unit	11.5	0.7	166	25	5	Positive
580	TG	TG(54:0) [NL-18:0]	FALSE	908.8	Unit	607.5	Unit	12.0	0.7	166	25	5	Positive
581	TG	TG(54:1) [NL-18:1]	FALSE	906.8	Unit	607.5	Unit	11.9	0.7	166	25	5	Positive
582	TG	TG(54:2) [NL-18:0]	FALSE	904.8	Unit	603.5	Unit	11.7	0.7	166	25	5	Positive
583	TG	TG(54:2) [NL-20:1]	FALSE	904.8	Unit	577.5	Unit	11.7	0.7	166	25	5	Positive
584	TG	TG(54:3) [NL-18:1]	FALSE	902.8	Unit	603.5	Unit	11.5	0.7	166	25	5	Positive
585	TG	TG(54:3) [NL-18:2]	FALSE	902.8	Unit	605.5	Unit	11.5	0.7	166	25	5	Positive
586	TG	TG(54:4) [NL-18:2]	FALSE	900.8	Unit	603.5	Unit	11.3	0.7	166	25	5	Positive
587	TG	TG(54:4) [NL-20:3]	FALSE	900.8	Unit	577.5	Unit	11.4	0.7	166	25	5	Positive
588	TG	TG(54:5) [NL-18:3]	FALSE	898.8	Unit	603.5	Unit	11.2	0.7	166	25	5	Positive
589	TG	TG(54:5) [NL-20:4]	FALSE	898.8	Unit	577.5	Unit	11.3	0.7	166	25	5	Positive
590	TG	TG(54:6) [NL-18:3]	FALSE	896.8	Unit	601.5	Unit	11.1	0.7	166	25	5	Positive
591	TG	TG(54:6) [NL-20:4]	FALSE	896.8	Unit	575.5	Unit	11.1	0.7	166	25	5	Positive
592	TG	TG(54:6) [NL-20:5]	FALSE	896.8	Unit	577.5	Unit	11.1	0.7	166	25	5	Positive
593	TG	TG(54:6) [NL-22:6]	FALSE	896.8	Unit	551.5	Unit	11.2	0.7	166	25	5	Positive
594	TG	TG(54:7) [NL-20:5]	FALSE	894.8	Unit	575.5	Unit	11.0	0.7	166	25	5	Positive
595	TG	TG(54:7) [NL-22:6]	FALSE	894.8	Unit	549.5	Unit	11.0	0.7	166	25	5	Positive
596	TG	TG(56:6) [NL-20:4]	FALSE	924.8	Unit	603.5	Unit	11.3	0.7	166	25	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	$\Delta RT$	碎裂电压	碰撞能量	池加速电压	极性
597	TG	TG(56:6) [NL-22:5]	FALSE	924.8	Unit	577.5	Unit	11.3	0.7	166	25	5	Positive
598	TG	TG(56:7) [NL-20:4]	FALSE	922.8	Unit	601.5	Unit	11.1	0.7	166	25	5	Positive
599	TG	TG(56:7) [NL-20:5]	FALSE	922.8	Unit	603.5	Unit	11.1	0.7	166	25	5	Positive
600	TG	TG(56:7) [NL-22:5]	FALSE	922.8	Unit	575.5	Unit	11.1	0.7	166	25	5	Positive
601	TG	TG(56:7) [NL-22:6]	FALSE	922.8	Unit	577.5	Unit	11.2	0.7	166	25	5	Positive
602	TG	TG(56:8) [NL-20:4]	FALSE	920.8	Unit	599.5	Unit	11.0	0.7	166	25	5	Positive
603	TG	TG(56:8) [NL-20:5]	FALSE	920.8	Unit	601.5	Unit	11.0	0.7	166	25	5	Positive
604	TG	TG(56:8) [NL-22:6]	FALSE	920.8	Unit	575.5	Unit	11.1	0.7	166	25	5	Positive
605	TG	TG(56:9) [NL-22:6]	FALSE	918.8	Unit	573.5	Unit	10.9	0.7	166	25	5	Positive
606	TG	TG(58:10) [NL-22:6]	FALSE	944.9	Unit	599.5	Unit	11.0	0.7	166	25	5	Positive
607	TG	TG(58:8) [NL-22:6]	FALSE	948.8	Unit	603.5	Unit	11.2	0.7	166	25	5	Positive
608	TG	TG(58:9) [NL-22:6]	FALSE	946.9	Unit	601.5	Unit	11.1	0.7	166	25	5	Positive
609	TG SIM	TG(48:0) [SIM]	FALSE	824.8	Unit	824.8	Unit	11.4	0.7	166	0	5	Positive
610	TG SIM	TG(48:1) [SIM]	FALSE	822.8	Unit	822.8	Unit	11.3	0.7	166	0	5	Positive
611	TG SIM	TG(48:2) [SIM]	FALSE	820.8	Unit	820.8	Unit	11.1	0.7	166	0	5	Positive
612	TG SIM	TG(48:3) [SIM]	FALSE	818.8	Unit	818.8	Unit	11.0	0.7	166	0	5	Positive
613	TG SIM	TG(49:1) [SIM]	FALSE	836.8	Unit	836.8	Unit	11.4	0.7	166	0	5	Positive
614	TG SIM	TG(50:0) [SIM]	FALSE	852.8	Unit	852.8	Unit	11.6	0.7	166	0	5	Positive
615	TG SIM	TG(50:1) [SIM]	FALSE	850.8	Unit	850.8	Unit	11.5	0.7	166	0	5	Positive
616	TG SIM	TG(50:2) [SIM]	FALSE	848.8	Unit	848.8	Unit	11.3	0.7	166	0	5	Positive
617	TG SIM	TG(50:3) [SIM]	FALSE	846.8	Unit	846.8	Unit	11.1	0.7	166	0	5	Positive
618	TG SIM	TG(50:4) [SIM]	FALSE	844.8	Unit	844.8	Unit	11.0	0.7	166	0	5	Positive
619	TG SIM	TG(51:0) [SIM]	FALSE	866.7	Unit	866.7	Unit	11.7	0.7	166	0	5	Positive
620	TG SIM	TG(51:1) [SIM]	FALSE	864.8	Unit	864.8	Unit	11.5	0.7	166	0	5	Positive
621	TG SIM	TG(51:2) [SIM]	FALSE	862.8	Unit	862.8	Unit	11.4	0.7	166	0	5	Positive
622	TG SIM	TG(52:1) [SIM]	FALSE	878.8	Unit	878.8	Unit	11.7	0.7	166	0	5	Positive
623	TG SIM	TG(52:2) [SIM]	FALSE	876.8	Unit	876.8	Unit	11.5	0.7	166	0	5	Positive
624	TG SIM	TG(52:3) [SIM]	FALSE	874.8	Unit	874.8	Unit	11.3	0.7	166	0	5	Positive
625	TG SIM	TG(52:4) [SIM]	FALSE	872.8	Unit	872.8	Unit	11.2	0.7	166	0	5	Positive
626	TG SIM	TG(52:5) [SIM]	FALSE	870.8	Unit	870.8	Unit	11.0	0.7	166	0	5	Positive
627	TG SIM	TG(53:2) [SIM]	FALSE	890.8	Unit	890.8	Unit	11.5	0.7	166	0	5	Positive
628	TG SIM	TG(54:0) [SIM]	FALSE	908.8	Unit	908.8	Unit	12.0	0.7	166	0	5	Positive
629	TG SIM	TG(54:1) [SIM]	FALSE	906.8	Unit	906.8	Unit	11.9	0.7	166	0	5	Positive
630	TG SIM	TG(54:2) [SIM]	FALSE	904.8	Unit	904.8	Unit	11.7	0.7	166	0	5	Positive
631	TG SIM	TG(54:3) [SIM]	FALSE	902.8	Unit	902.8	Unit	11.5	0.7	166	0	5	Positive
632	TG SIM	TG(54:4) [SIM]	FALSE	900.8	Unit	900.8	Unit	11.3	0.7	166	0	5	Positive
633	TG SIM	TG(54:5) [SIM]	FALSE	898.8	Unit	898.8	Unit	11.2	0.7	166	0	5	Positive
634	TG SIM	TG(54:6) [SIM]	FALSE	896.8	Unit	896.8	Unit	11.1	0.7	166	0	5	Positive
635	TG SIM	TG(54:7) [SIM]	FALSE	894.8	Unit	894.8	Unit	11.0	0.7	166	0	5	Positive
636	TG SIM	TG(56:6) [SIM]	FALSE	924.8	Unit	924.8	Unit	11.3	0.7	166	0	5	Positive
637	TG SIM	TG(56:7) [SIM]	FALSE	922.8	Unit	922.8	Unit	11.2	0.7	166	0	5	Positive
638	TG SIM	TG(56:8) [SIM]	FALSE	920.8	Unit	920.8	Unit	11.1	0.7	166	0	5	Positive
639	TG SIM	TG(56:9) [SIM]	FALSE	918.8	Unit	918.8	Unit	10.9	0.7	166	0	5	Positive
640	TG SIM	TG(58:10) [SIM]	FALSE	944.9	Unit	944.9	Unit	11.0	0.7	166	0	5	Positive
641	TG SIM	TG(O-50:1) [SIM]	FALSE	836.8	Unit	836.8	Unit	11.8	0.7	166	0	5	Positive
642	TG SIM	TG(O-50:2) [SIM]	FALSE	834.8	Unit	834.8	Unit	11.6	0.7	166	0	5	Positive

化合物编号	化合物组	化合物名称	是否为内标?	母离子	MS1分辨率	子离子	MS2分辨率	保留时间 (min)	ΔRT	碎裂电压	碰撞能量	池加速电压	极性
643	TG SIM	TG(O-52:1) [SIM]	FALSE	864.8	Unit	864.8	Unit	12.0	0.7	166	0	5	Positive
644	TG SIM	TG(O-52:2) [SIM]	FALSE	862.8	Unit	862.8	Unit	11.8	0.7	166	0	5	Positive
645	TG SIM	TG(O-54:2) [SIM]	FALSE	890.8	Unit	890.8	Unit	12.0	0.7	166	0	5	Positive
646	TG (O)	TG(O-50:1) [NL-15:0]	FALSE	836.8	Unit	577.5	Unit	11.8	0.7	166	25	5	Positive
647	TG (O)	TG(O-50:1) [NL-16:0]	FALSE	836.8	Unit	563.5	Unit	11.8	0.7	166	25	5	Positive
648	TG (O)	TG(O-50:1) [NL-17:1]	FALSE	836.8	Unit	551.5	Unit	11.8	0.7	166	25	5	Positive
649	TG (O)	TG(O-50:1) [NL-18:1]	FALSE	836.8	Unit	537.5	Unit	11.8	0.7	166	25	5	Positive
650	TG (O)	TG(O-50:2) [NL-16:1]	FALSE	834.8	Unit	563.5	Unit	11.6	0.7	166	25	5	Positive
651	TG (O)	TG(O-50:2) [NL-18:1]	FALSE	834.8	Unit	535.5	Unit	11.6	0.7	166	25	5	Positive
652	TG (O)	TG(O-50:2) [NL-18:2]	FALSE	834.8	Unit	537.5	Unit	11.6	0.7	166	25	5	Positive
653	TG (O)	TG(O-52:0) [NL-16:0]	FALSE	866.8	Unit	593.5	Unit	12.2	0.7	166	25	5	Positive
654	TG (O)	TG(O-52:1) [NL-16:0]	FALSE	864.8	Unit	591.5	Unit	12.0	0.7	166	25	5	Positive
655	TG (O)	TG(O-52:1) [NL-18:1]	FALSE	864.8	Unit	565.5	Unit	12.0	0.7	166	25	5	Positive
656	TG (O)	TG(O-52:2) [NL-16:0]	FALSE	862.8	Unit	589.5	Unit	11.8	0.7	166	25	5	Positive
657	TG (O)	TG(O-52:2) [NL-17:1]	FALSE	862.8	Unit	577.5	Unit	11.8	0.7	166	25	5	Positive
658	TG (O)	TG(O-52:2) [NL-18:1]	FALSE	862.8	Unit	563.5	Unit	11.8	0.7	166	25	5	Positive
659	TG (O)	TG(O-54:2) [NL-17:1]	FALSE	890.8	Unit	605.5	Unit	12.0	0.7	166	25	5	Positive
660	TG (O)	TG(O-54:2) [NL-18:1]	FALSE	890.8	Unit	591.5	Unit	12.0	0.7	166	25	5	Positive
661	TG (O)	TG(O-54:3) [NL-17:1]	FALSE	888.8	Unit	603.5	Unit	11.8	0.7	166	25	5	Positive
662	TG (O)	TG(O-54:3) [NL-18:1]	FALSE	888.8	Unit	589.5	Unit	11.8	0.7	166	25	5	Positive
663	TG (O)	TG(O-54:4) [NL-17:1]	FALSE	886.8	Unit	601.5	Unit	11.6	0.7	166	25	5	Positive
664	TG (O)	TG(O-54:4) [NL-18:2]	FALSE	886.8	Unit	589.5	Unit	11.6	0.7	166	25	5	Positive
665	Ubiquinone	Ubiquinone	FALSE	880.7	Unit	197	Unit	11.0	0.4	166	17	5	Positive

<sup>1</sup> 此处的离子对数量少于脂质种类的最终列表, 因为某些离子对可以监测在积分时分离的多个峰

表 A2. 每种脂质的推荐参比化合物以及响应因子和 %CV 值

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
AC(10:0)	316.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	7.2	<1
AC(12:0)	344.3 → 85.1	AC(12:0) [reference]	AC(16:0) d3	1	2.7	<1
AC(12:1)	342.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	5.4	<1
AC(13:0)	358.3 → 85.1	AC(12:0) [reference]	AC(16:0) d3	1	1.8	1.16
AC(14:0)	372.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	1.7	2.02
AC(14:0)-OH	388.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	4.0	1.38
AC(14:1)	370.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	2.6	<1
AC(14:1)-OH	386.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	5.9	<1
AC(14:2)	368.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	3.1	<1
AC(15:0) (a)	386.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	6.2	2.06
AC(15:0) (b)	386.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	10.3	12.37
AC(16:0)	400.4 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	1.6	<1
AC(16:0)-OH	416.4 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	7.9	3.41
AC(16:1)	398.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	2.5	1.29
AC(16:1)-OH	414.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	6.1	6.41
AC(17:0) (a)	414.4 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	6.7	7.62
AC(17:0) (b)	414.4 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	6.2	8.87

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
AC(18:0)	428.4 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	2.8	<1
AC(18:0)-OH	444.4 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	11.9	8.3
AC(18:1)	426.4 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	1.9	<1
AC(18:1)-OH	442.4 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	7.9	1.29
AC(18:2)	424.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	2.4	<1
AC(18:3)	422.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	4.6	7.71
AC(20:3) (a)	450.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	8.0	6.11
AC(20:3) (b)	450.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	17.1	23.07
AC(20:3)-OH	466.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	11.4	<1
AC(20:4)	448.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	14.1	7.47
AC(20:5)	446.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	8.2	2.97
AC(22:5)	474.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	17.9	<1
AC(22:5)-OH	490.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	7.9	1.24
AC(22:6)	472.3 → 85.1	AcylCarnitine 16:0 d3 (IS) [reference]	AC(16:0) d3	1	16.8	24.67
AC(24:0)	512.3 → 85.1	AC(24:0) [reference]	AC(16:0) d3	1	3.0	<1
AC(24:1) (a)	510.3 → 85.1	AC(24:0) [reference]	AC(16:0) d3	1	5.9	2.21
AC(24:1) (b)	510.3 → 85.1	AC(24:0) [reference]	AC(16:0) d3	1	9.8	6.96
AC(26:0)	540.3 → 85.1	AC(24:0) [reference]	AC(16:0) d3	1	3.3	<1
AC(26:1)	538.3 → 85.1	AC(24:0) [reference]	AC(16:0) d3	1	3.7	<1
CA	426.3 → 355.3	Cholic Acid d4 (IS) [reference]	Cholic acid d4	1	5.4	10
CE(14:0)	614.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	4.3	18.4	<1
CE(15:0)	628.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	2.4	12.0	1.03
CE(16:0)	642.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1.7	2.3	<1
CE(16:1)	640.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1	10.4	<1
CE(16:2)	638.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.5	9.8	<1
CE(17:0)	656.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1.3	4.9	<1
CE(17:1)	654.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.7	7.5	<1
CE(18:0)	670.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1.1	2.0	<1
CE(18:1)	668.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.6	3.1	<1
CE(18:2)	666.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.3	4.5	<1
CE(18:2) [+OH]	682.6 → 369.3	CE(18:3) [reference]	CE(18:0) d6	1	9.4	<1
CE(18:3)	664.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.3	7.9	<1
CE(20:0)	698.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1.1	12.9	6.67
CE(20:1)	696.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.4	5.1	3.45
CE(20:2)	694.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.2	2.8	<1
CE(20:3)	692.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.2	2.9	<1
CE(20:4)	690.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.2	4.1	<1
CE(20:4) [+OH]	706.6 → 369.3	CE(20:5) [reference]	CE(18:0) d6	1	10.9	<1
CE(20:5)	688.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.2	3.7	<1
CE(22:0)	726.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.6	14.8	8.7
CE(22:1)	724.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.3	17.0	3.71
CE(22:4)	718.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.2	4.4	<1
CE(22:5)	716.6 → 369.3	CE(22:6) [reference]	CE(18:0) d6	0.2	4.3	<1
CE(22:6)	714.6 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.2	6.1	<1
CE(24:0)	754.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.5	18.7	16.71
CE(24:1)	752.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.3	16.6	6.78
CE(24:4)	746.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.1	8.4	<1

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
CE(24:5)	744.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.1	8.1	<1
CE(24:6)	742.7 → 369.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	0.1	13.7	<1
Cer(d16:1/16:0)	510.6 → 236.3	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	12.6	1.31
Cer(d16:1/18:0)	538.6 → 236.3	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	7.5	3.53
Cer(d16:1/20:0)	566.6 → 236.3	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	4.2	2.12
Cer(d16:1/22:0)	594.6 → 236.3	Cer(d16:1/22:0) [reference]	Cer(d18:1-d7/18:0)	1	14.3	1.33
Cer(d16:1/23:0)	608.6 → 236.3	Cer(d16:1/23:0) [reference]	Cer(d18:1-d7/18:0)	1	14.1	1.7
Cer(d16:1/24:0)	622.6 → 236.3	Cer(d16:1/24:0) [reference]	Cer(d18:1-d7/18:0)	1	15.3	1.2
Cer(d16:1/24:1)	620.6 → 236.3	Cer(d16:1/24:1) [reference]	Cer(d18:1-d7/18:0)	1	14.2	<1
Cer(d17:1/16:0)	524.6 → 250.3	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	12.9	7.44
Cer(d17:1/18:0)	552.6 → 250.3	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	11.1	28.25
Cer(d17:1/20:0)	580.6 → 250.3	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	8.4	8.01
Cer(d17:1/22:0)	608.6 → 250.3	Cer(d17:1/22:0) [reference]	Cer(d18:1-d7/18:0)	1	13.5	2.66
Cer(d17:1/23:0)	622.6 → 250.3	Cer(d17:1/23:0) [reference]	Cer(d18:1-d7/18:0)	1	16.2	2.04
Cer(d17:1/24:0)	636.6 → 250.3	Cer(d17:1/24:0) [reference]	Cer(d18:1-d7/18:0)	1	14.5	2.13
Cer(d17:1/24:1)	634.6 → 250.3	Cer(d17:1/24:1) [reference]	Cer(d18:1-d7/18:0)	1	14.2	<1
Cer(d18:1/16:0)	538.5 → 264.3	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	3.9	<1
Cer(d18:1/18:0)	566.6 → 264.3	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	3.1	1.11
Cer(d18:1/20:0)	594.6 → 264.3	Cer(d18:1/21:0) [reference]	Cer(d18:1-d7/18:0)	1	13.7	3.03
Cer(d18:1/21:0)	608.6 → 264.3	Cer(d18:1/21:0) [reference]	Cer(d18:1-d7/18:0)	1	14.2	2.95
Cer(d18:1/22:0)	622.6 → 264.3	Cer(d18:1/22:0) [reference]	Cer(d18:1-d7/18:0)	1	13.5	<1
Cer(d18:1/23:0)	636.6 → 264.3	Cer(d18:1/23:0) [reference]	Cer(d18:1-d7/18:0)	1	13.4	<1
Cer(d18:1/24:0)	650.6 → 264.3	Cer(d18:1/24:0) [reference]	Cer(d18:1-d7/18:0)	1	15.9	<1
Cer(d18:1/24:1)	648.6 → 264.3	Cer(d18:1/24:1) [reference]	Cer(d18:1-d7/18:0)	1	13.5	<1
Cer(d18:1/26:0)	678.6 → 264.3	Cer(d18:1/26:0) [reference]	Cer(d18:1-d7/18:0)	1	17.3	13.36
Cer(d18:2/16:0)	536.5 → 262.3	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	9.9	1.09
Cer(d18:2/18:0)	564.6 → 262.3	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	7.2	<1
Cer(d18:2/20:0)	592.6 → 262.3	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	7.4	<1
Cer(d18:2/22:0)	620.6 → 262.3	Cer(d18:2/22:0) [reference]	Cer(d18:1-d7/18:0)	1	14.5	<1
Cer(d18:2/23:0)	634.6 → 262.3	Cer(d18:2/23:0) [reference]	Cer(d18:1-d7/18:0)	1	14.4	<1
Cer(d18:2/24:0)	648.6 → 262.3	Cer(d18:2/24:0) [reference]	Cer(d18:1-d7/18:0)	1	16.0	<1
Cer(d18:2/24:1)	646.6 → 262.3	Cer(d18:2/24:1) [reference]	Cer(d18:1-d7/18:0)	1	14.4	<1
Cer(d18:2/26:0)	676.6 → 262.3	Cer(d18:2/24:0) [reference]	Cer(d18:1-d7/18:0)	1	18.2	2.79
Cer(d19:1/18:0)	580.6 → 278.3	Cer(d19:1/20:0) [reference]	Cer(d18:1-d7/18:0)	1	13.5	4.15
Cer(d19:1/20:0)	608.6 → 278.3	Cer(d19:1/20:0) [reference]	Cer(d18:1-d7/18:0)	1	16.3	13.82
Cer(d19:1/22:0)	636.6 → 278.3	Cer(d19:1/22:0) [reference]	Cer(d18:1-d7/18:0)	1	15.2	<1
Cer(d19:1/23:0)	650.6 → 278.3	Cer(d19:1/23:0) [reference]	Cer(d18:1-d7/18:0)	1	15.4	<1
Cer(d19:1/24:0)	664.6 → 278.3	Cer(d19:1/24:0) [reference]	Cer(d18:1-d7/18:0)	1	16.0	<1
Cer(d19:1/24:1)	662.6 → 278.3	Cer(d19:1/24:1) [reference]	Cer(d18:1-d7/18:0)	1	14.8	<1
Cer(d20:1/22:0)	650.6 → 292.3	Cer(d20:1/22:0) [reference]	Cer(d18:1-d7/18:0)	1	17.1	8.75
Cer(d20:1/23:0)	664.6 → 292.3	Cer(d20:1/23:0) [reference]	Cer(d18:1-d7/18:0)	1	17.0	5.41
Cer(d20:1/24:0)	678.6 → 292.3	Cer(d20:1/24:0) [reference]	Cer(d18:1-d7/18:0)	1	16.6	5.37
Cer(d20:1/24:1)	676.6 → 292.3	Cer(d20:1/24:1) [reference]	Cer(d18:1-d7/18:0)	1	16.4	1.26
Cer(m18:0/20:0)	580.6 → 268.4	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	14.2	3
Cer(m18:0/22:0)	608.6 → 268.4	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	15.3	<1
Cer(m18:0/23:0)	622.6 → 268.4	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	14.8	1.08
Cer(m18:0/24:0)	636.6 → 268.4	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	15.7	<1

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
Cer(m18:0/24:1)	634.6 → 268.4	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	15.2	1.36
Cer(m18:1/18:0)	550.6 → 266.4	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	14.2	<1
Cer(m18:1/20:0)	578.6 → 266.4	Cer(m18:1/20:0) [reference]	Cer(d18:1-d7/18:0)	1	14.6	<1
Cer(m18:1/22:0)	606.6 → 266.4	Cer(m18:1/20:0) [reference]	Cer(d18:1-d7/18:0)	1	14.4	<1
Cer(m18:1/23:0)	620.6 → 266.4	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	13.2	<1
Cer(m18:1/24:0)	634.6 → 266.4	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	15.3	<1
Cer(m18:1/24:1)	632.6 → 266.4	Cer(d18:1-d7/18:0) (IS) [reference]	Cer(d18:1-d7/18:0)	1	14.4	<1
Cer1P(d18:1/16:0)	618.4 → 264.3	Cer1P(d18:1/16:0) [reference]	Cer(d18:1-d7/18:0)	1	16.6	<1
COH	369.4 → 161.2	COH-d7 (IS) [reference]	COH(d7)	1	7.9	<1
DE(16:0)	640.6 → 367.4	DE(18:1) ester d6 (IS) [reference]	CE(18:0) d6	1	16.7	1.74
DE(18:1)	666.6 → 367.4	DE(18:1) ester d6 (IS) [reference]	CE(18:0) d6	1	6.8	<1
DE(18:2)	664.6 → 367.4	DE(18:1) ester d6 (IS) [reference]	CE(18:0) d6	1	7.7	<1
DE(20:4)	688.6 → 367.4	DE(18:1) ester d6 (IS) [reference]	CE(18:0) d6	1	7.8	<1
DG(14:0_16:0)	558.5 → 313.3	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	3.3	28.6
DG(14:0_18:2)	582.5 → 285.2	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	7.8	1.94
DG(16:0_16:0)	586.5 → 313.2	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	0.5	4.5	16.01
DG(16:0_16:1)	584.5 → 313.2	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	2.9	10.96
DG(16:0_18:1)	612.6 → 313.3	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	4.2	2.83
DG(16:0_18:2)	610.5 → 313.2	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	4.8	<1
DG(16:0_20:4)	634.5 → 313.2	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	4.1	6.05
DG(16:0_22:5)	660.6 → 313.3	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	7.7	7.88
DG(16:0_22:6)	658.5 → 313.2	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	7.2	4.45
DG(16:1_18:1)	610.5 → 339.2	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	2.7	3.86
DG(18:0_18:1)	640.6 → 341.3	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	3.6	5.13
DG(18:0_18:2)	638.6 → 341.3	DG(18:0_18:2) [reference]	DG(15:0_18:1) d7	1	4.3	1.73
DG(18:0_20:4)	662.6 → 341.3	DG(18:0_20:4) [reference]	DG(15:0_18:1) d7	1	8.7	1.64
DG(18:1_18:1)	638.6 → 339.3	DG(18:0_18:2) [reference]	DG(15:0_18:1) d7	1	3.4	1.2
DG(18:1_18:2)	636.6 → 339.3	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	2.4	<1
DG(18:1_18:3)	634.5 → 339.2	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	4.0	<1
DG(18:1_20:3)	662.6 → 339.3	DG(18:0_20:4) [reference]	DG(15:0_18:1) d7	1	4.7	1.22
DG(18:1_20:4)	660.6 → 339.3	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	2.6	<1
DG(18:1_20:5)	658.6 → 339.3	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	3.5	2.71
DG(18:1_22:5)	686.6 → 339.3	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	5.1	3.15
DG(18:1_22:6)	684.6 → 339.3	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	3.6	1.27
DG(18:2_18:2)	634.5 → 337.2	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	0.5	3.0	<1
DG(18:2_20:4)	658.5 → 337.2	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	3.9	<1
DG(18:2_22:6)	682.6 → 337.3	DG(15:0 18:1) d7 (IS) [reference]	DG(15:0_18:1) d7	1	4.8	<1
dhCer(d18:0/16:0)	540.5 → 284.3	dhCer(d18:0/8:0) [reference]	dhCer(d18:0/8:0)	1	11.8	32.71
dhCer(d18:0/18:0)	568.6 → 284.3	dhCer(d18:0/8:0) [reference]	dhCer(d18:0/8:0)	1	9.4	23.82
dhCer(d18:0/20:0)	596.6 → 284.3	dhCer(d18:0/8:0) [reference]	dhCer(d18:0/8:0)	1	14.6	16.64
dhCer(d18:0/22:0)	624.6 → 284.3	dhCer(d18:0/8:0) [reference]	dhCer(d18:0/8:0)	1	15.4	1.67
dhCer(d18:0/24:0)	652.7 → 284.3	dhCer(d18:0/8:0) [reference]	dhCer(d18:0/8:0)	1	16.1	2.87
dhCer(d18:0/24:1)	650.6 → 284.3	dhCer(d18:0/8:0) [reference]	dhCer(d18:0/8:0)	1	14.3	<1
dimethyl-CE(18:1)	696.6 → 397.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1	6.0	<1
dimethyl-CE(18:2)	694.6 → 397.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1	2.6	<1
dimethyl-CE(20:4)	718.6 → 397.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1	5.0	<1
dxCA	410.3 → 357.3	Cholic Acid d4 (IS) [reference]	Cholic acid d4	1	4.7	8.07



化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
FA(14:0)	227.2 → 227.2	FA(22:6) d5 (IS) [reference]	FFA(18:1) d9	1	6.2	67.45
FA(16:0)	255.2 → 255.2	FA(18:1) d9 (IS) [reference]	FFA(18:1) d9	1	2.0	23.4
FA(16:1)	253.2 → 253.2	FA(20:4) d11 (IS) [reference]	FFA(18:1) d9	1	4.8	14.87
FA(17:0)	269.2 → 269.2	FA(18:1) d9 (IS) [reference]	FFA(18:1) d9	1	3.6	79.9
FA(17:1)	267.2 → 267.2	FA(18:1) d9 (IS) [reference]	FFA(18:1) d9	1	2.8	70.54
FA(18:0)	283.3 → 283.3	FA(18:1) d9 (IS) [reference]	FFA(18:1) d9	1	4.5	94.5
FA(18:1)	281.2 → 281.2	FA(18:1) d9 (IS) [reference]	FFA(18:1) d9	1	1.6	2.23
FA(18:2)	279.2 → 279.2	FA(18:1) d9 (IS) [reference]	FFA(18:1) d9	1	3.1	10.48
FA(18:3)	277.2 → 277.2	FA(22:6) d5 (IS) [reference]	FFA(18:1) d9	1	3.8	9.42
FA(20:2)	307.3 → 307.3	FA(18:1) d9 (IS) [reference]	FFA(18:1) d9	1	6.7	73.72
FA(20:3)	305.2 → 305.2	FA(18:1) d9 (IS) [reference]	FFA(18:1) d9	1	3.9	76.41
FA(20:4)	303.2 → 303.2	FA(20:4) d11 (IS) [reference]	FFA(18:1) d9	1	5.1	14.41
FA(20:5)	301.2 → 301.2	FA(22:6) d5 (IS) [reference]	FFA(18:1) d9	1	7.3	26.16
FA(22:4)	331.3 → 331.3	FA(18:1) d9 (IS) [reference]	FFA(18:1) d9	1	4.5	34.32
FA(22:5)	329.2 → 329.2	FA(20:4) d11 (IS) [reference]	FFA(18:1) d9	1	7.8	43.67
FA(22:6)	327.2 → 327.2	FA(22:6) d5 (IS) [reference]	FFA(18:1) d9	1	4.6	26.58
GM3(d18:1/16:0)	1153.7 → 264.3	GM3(d18:1/16:0) [reference]	Hex3Cer(d18:1/17:0)	1	4.2	<1
GM3(d18:1/18:0)	1181.8 → 264.3	GM3(d18:1/16:0) [reference]	Hex3Cer(d18:1/17:0)	1	4.4	<1
GM3(d18:1/20:0)	1209.8 → 264.3	GM3(d18:1/16:0) [reference]	Hex3Cer(d18:1/17:0)	1	7.4	<1
GM3(d18:1/22:0)	1237.8 → 264.3	GM3(d18:1/22:0) [reference]	Hex3Cer(d18:1/17:0)	1	5.6	<1
GM3(d18:1/24:0)	1265.8 → 264.3	GM3(d18:1/24:0) [reference]	Hex3Cer(d18:1/17:0)	1	4.9	<1
GM3(d18:1/24:1)	1263.8 → 264.3	GM3(d18:1/22:0) [reference]	Hex3Cer(d18:1/17:0)	1	5.1	<1
Hex2Cer(d16:1/16:0)	834.6 → 236.3	Hex2Cer(d18:1/15:0) d7 (IS) [reference]	Hex2Cer(d18:1/15:0) d7	1	4.2	<1
Hex2Cer(d16:1/24:1)	944.7 → 236.3	Hex2Cer(d16:1/24:1) [reference]	Hex2Cer(d18:1/15:0) d7	1	8.6	<1
Hex2Cer(d18:1/16:0)	862.6 → 264.3	Hex2Cer(d18:1/16:0) [reference]	Hex2Cer(d18:1/15:0) d7	1	2.4	<1
Hex2Cer(d18:1/20:0)	918.7 → 264.3	Hex2Cer(d18:1/22:0) [reference]	Hex2Cer(d18:1/15:0) d7	1	9.0	<1
Hex2Cer(d18:1/22:0)	946.7 → 264.3	Hex2Cer(d18:1/22:0) [reference]	Hex2Cer(d18:1/15:0) d7	1	3.8	<1
Hex2Cer(d18:1/24:0)	974.8 → 264.3	Hex2Cer(d18:1/24:0) [reference]	Hex2Cer(d18:1/15:0) d7	1	4.4	<1
Hex2Cer(d18:1/24:1)	972.7 → 264.3	Hex2Cer(d18:1/24:1) [reference]	Hex2Cer(d18:1/15:0) d7	1	2.6	<1
Hex2Cer(d18:2/16:0)	860.6 → 262.3	Hex2Cer(d18:1/15:0) d7 (IS) [reference]	Hex2Cer(d18:1/15:0) d7	1	4.6	<1
Hex2Cer(d18:2/24:1)	970.7 → 262.3	Hex2Cer(d18:2/24:1) [reference]	Hex2Cer(d18:1/15:0) d7	1	4.7	<1
Hex3Cer(d18:1/16:0)	1024.7 → 264.3	Hex3Cer(d18:1/17:0) (IS) [reference]	Hex3Cer(d18:1/17:0)	1	4.4	<1
Hex3Cer(d18:1/18:0)	1052.7 → 264.3	Hex3Cer(d18:1/17:0) (IS) [reference]	Hex3Cer(d18:1/17:0)	1	11.2	1.58
Hex3Cer(d18:1/22:0)	1108.8 → 264.3	Hex3Cer(d18:1/17:0) (IS) [reference]	Hex3Cer(d18:1/17:0)	1	5.3	<1
Hex3Cer(d18:1/24:0)	1136.8 → 264.3	Hex3Cer(d18:1/17:0) (IS) [reference]	Hex3Cer(d18:1/17:0)	1	5.3	<1
Hex3Cer(d18:1/24:1)	1134.8 → 264.3	Hex3Cer(d18:1/17:0) (IS) [reference]	Hex3Cer(d18:1/17:0)	1	5.1	<1
HexCer(d16:1/18:0)	700.6 → 236.3	HexCer(d16:1/20:0) [reference]	HexCer(d18:1/15:0) d7	1	16.7	3.78
HexCer(d16:1/20:0)	728.6 → 236.3	HexCer(d16:1/20:0) [reference]	HexCer(d18:1/15:0) d7	1	9.8	2.88
HexCer(d16:1/22:0)	756.7 → 236.3	HexCer(d16:1/24:0) [reference]	HexCer(d18:1/15:0) d7	1	6.8	<1
HexCer(d16:1/24:0)	784.7 → 236.3	HexCer(d16:1/24:0) [reference]	HexCer(d18:1/15:0) d7	1	6.2	<1
HexCer(d18:1/16:0)	700.6 → 264.3	HexCer(d18:1/16:0) [reference]	HexCer(d18:1/15:0) d7	1	8.6	<1
HexCer(d18:1/18:0)	728.6 → 264.3	HexCer(d18:1/18:0) [reference]	HexCer(d18:1/15:0) d7	1	6.9	<1
HexCer(d18:1/20:0)	756.6 → 264.3	HexCer(d18:1/20:0) [reference]	HexCer(d18:1/15:0) d7	1	7.2	<1
HexCer(d18:1/22:0)	784.7 → 264.3	HexCer(d18:1/22:0) [reference]	HexCer(d18:1/15:0) d7	1	2.5	<1
HexCer(d18:1/24:0)	812.7 → 264.3	HexCer(d18:1/24:0) [reference]	HexCer(d18:1/15:0) d7	1	3.5	<1
HexCer(d18:1/24:1)	810.7 → 264.3	HexCer(d18:1/24:1) [reference]	HexCer(d18:1/15:0) d7	1	2.6	<1
HexCer(d18:2/20:0)	754.6 → 262.3	HexCer(d18:2/20:0) [reference]	HexCer(d18:1/15:0) d7	1	13.0	1.36

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
HexCer(d18:2/22:0)	782.7 → 262.3	HexCer(d18:2/20:0) [reference]	HexCer(d18:1/15:0) d7	1	7.1	<1
HexCer(d18:2/24:0)	810.7 → 262.3	HexCer(d18:2/24:0) [reference]	HexCer(d18:1/15:0) d7	1	4.2	<1
LPC(14:0) [sn1]	468.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	2.0	<1
LPC(14:0) [sn2]	468.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	3.5	<1
LPC(15:0) [sn1]	482.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	1.9	17.68
LPC(15:0) [sn2]	482.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	3.7	<1
LPC(15-MHDA) [sn1] / LPC(17:0) [sn2]	510.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	2.1	<1
LPC(15-MHDA) [sn1] [104_sn1]	510.4 → 104.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	3.0	<1
LPC(15-MHDA) [sn2]	510.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	3.7	<1
LPC(16:0) [sn1]	496.3 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	1.8	<1
LPC(16:0) [sn2]	496.3 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	2.1	<1
LPC(16:1) [sn1]	494.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	1.5	<1
LPC(16:1) [sn2]	494.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	4.4	<1
LPC(17:0) [sn1]	510.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	2.1	<1
LPC(17:1) (a) [sn1] [104_sn1]	508.4 → 104.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	3.3	<1
LPC(17:1) [sn1] (a) / LPC(17:1) [sn2] (b)	508.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	2.3	<1
LPC(17:1) [sn1] (b)	508.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	4.8	2.07
LPC(17:1) [sn2] (a)	508.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	5.6	<1
LPC(18:0) [sn1]	524.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	2.0	<1
LPC(18:0) [sn2]	524.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	2.8	<1
LPC(18:1) [sn1]	522.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	1.5	<1
LPC(18:1) [sn2]	522.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	1.9	<1
LPC(18:2) [+OH]	536.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	8.8	4.8
LPC(18:2) [sn1]	520.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	1.7	<1
LPC(18:2) [sn2]	520.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	3.6	<1
LPC(18:3) (a) [sn1] [104_sn1]	518.3 → 104.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	4.8	<1
LPC(18:3) [sn1] (a)/LPC(18:3) [sn2] (b)	518.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	3.4	<1
LPC(18:3) [sn1] (b)	518.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	3.7	<1
LPC(18:3) [sn2] (a)	518.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	5.3	<1
LPC(19:0) (a) [sn1] [104_sn1]	538.4 → 104.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	5.9	<1
LPC(19:0) [sn1] (a) / LPC(19:0) [sn2] (b)	538.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	4.3	1.21
LPC(19:0) [sn1] (b)	538.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	3.9	<1
LPC(19:0) [sn2] (a)	538.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	6.9	6.16
LPC(19:1) (a)	536.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	10.0	4.92
LPC(19:1) (b)	536.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	4.4	2.96
LPC(19:1) (c)	536.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	9.0	28.49
LPC(20:0) [sn1]	552.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	4.5	<1
LPC(20:0) [sn2]	552.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	4.8	2.53
LPC(20:1) [sn1]	550.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	4.1	<1
LPC(20:1) [sn2]	550.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	4.1	1
LPC(20:2) [sn1]	548.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	2.8	<1
LPC(20:2) [sn2]	548.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	4.1	<1
LPC(20:3) [sn1]	546.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	2.4	<1
LPC(20:3) [sn2]	546.4 → 184.1	LPC(0-16:0) [reference]	LPC(18:1) d7	1	3.3	<1
LPC(20:4) [+OH]	560.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	15.4	25.83
LPC(20:4) [sn1]	544.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	3.3	<1
LPC(20:4) [sn2]	544.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	4.0	<1

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
LPC(20:5) [sn1]	542.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	4.3	<1
LPC(20:5) [sn2]	542.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	5.6	<1
LPC(22:0) [sn1]	580.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	8.0	<1
LPC(22:0) [sn2]	580.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	8.6	<1
LPC(22:1) [sn1]	578.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	5.8	3.53
LPC(22:1) [sn2]	578.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	8.3	15.86
LPC(22:4) [sn1]	572.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	4.7	<1
LPC(22:4) [sn2]	572.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	5.5	<1
LPC(22:5) (n3) [sn1] [104_sn1]	570.4 → 104.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	3.3	<1
LPC(22:5) [sn1] (n3)/LPC(22:5) [sn2] (n6)	570.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	3.0	<1
LPC(22:5) [sn1] (n6)	570.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	4.8	<1
LPC(22:5) [sn2] (n3)	570.4 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	4.0	<1
LPC(22:6) [sn1]	568.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	3.8	<1
LPC(22:6) [sn2]	568.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	5.0	<1
LPC(22:6) [+OH]	584.3 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	38.1	53.47
LPC(24:0) [sn1]	608.5 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	6.9	<1
LPC(24:0) [sn2]	608.5 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	8.0	<1
LPC(26:0) [sn1]	636.5 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	6.4	43.83
LPC(26:0) [sn2]	636.5 → 184.1	LPC(18:1) d7 (IS) [reference]	LPC(18:1) d7	1	7.3	51.49
LPC(O-16:0)	482.4 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	2.1	<1
LPC(O-18:0)	510.4 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	3.5	<1
LPC(O-18:1)	508.4 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	2.4	<1
LPC(O-20:0)	538.4 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	6.3	<1
LPC(O-20:1)	536.4 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	4.3	<1
LPC(O-22:0)	566.5 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	5.8	<1
LPC(O-22:1)	564.4 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	7.1	<1
LPC(O-24:0)	594.5 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	4.4	<1
LPC(O-24:1)	592.5 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	5.6	1.54
LPC(O-24:2)	590.5 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	7.9	<1
LPC(P-16:0)	480.3 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	2.2	<1
LPC(P-17:0) (a)	494.3 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	5.3	3.84
LPC(P-17:0) (b)	494.3 → 104.1	LPC(O-16:0) [reference]	LPC(18:1) d7	1	6.7	8.79
LPC(P-18:0)	508.3 → 104.1	LPC(P-18:0) [reference]	LPC(18:1) d7	1	5.1	<1
LPC(P-18:1)	506.3 → 104.1	LPC(P-18:1) [reference]	LPC(18:1) d7	1	4.6	<1
LPC(P-20:0)	536.3 → 104.1	LPC(P-18:1) [reference]	LPC(18:1) d7	1	11.3	<1
LPE(16:0) [sn1]	454.3 → 313.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	5.5	<1
LPE(16:0) [sn2]	454.3 → 313.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	5.1	<1
LPE(17:0)	468.3 → 327.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	22.8	7.3
LPE(18:0) [sn1]	482.3 → 341.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	3.6	<1
LPE(18:0) [sn2]	482.3 → 341.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	5.3	<1
LPE(18:1) [sn1]	480.3 → 339.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	5.0	<1
LPE(18:1) [sn2]	480.3 → 339.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	5.6	<1
LPE(18:2) [sn1]	478.3 → 337.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	3.9	<1
LPE(18:2) [sn2]	478.3 → 337.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	5.0	<1
LPE(20:4) [sn1]	502.3 → 361.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	5.3	<1
LPE(20:4) [sn2]	502.3 → 361.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	6.1	<1
LPE(22:6) [sn1]	526.3 → 385.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	5.0	<1

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
LPE(22:6) [sn2]	526.3 → 385.3	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	7.3	<1
LPE(P-16:0)	438.3 → 266.4	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	11.2	<1
LPE(P-18:0)	466.3 → 294.4	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	9.6	<1
LPE(P-18:1)	464.3 → 292.4	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	15.8	<1
LPE(P-20:0)	494.3 → 322.4	LPE(18:1) d7 (IS) [reference]	LPE(18:1) d7	1	21.6	21.11
LPI(18:0)	618.3 → 341.3	LPI(18:0) [reference]	LPI(13:0)	1	18.7	2.56
LPI(18:1) [sn1]	616.3 → 339.3	LPI 13:0 (IS) [reference]	LPI(13:0)	1	10.8	<1
LPI(18:1) [sn2]	616.3 → 339.3	LPI 13:0 (IS) [reference]	LPI(13:0)	1	25.4	<1
LPI(18:2) [sn1]	614.3 → 337.3	LPI 13:0 (IS) [reference]	LPI(13:0)	1	10.6	<1
LPI(18:2) [sn2]	614.3 → 337.3	LPI 13:0 (IS) [reference]	LPI(13:0)	1	10.6	3.74
LPI(20:4) [sn1]	638.3 → 361.3	LPI 13:0 (IS) [reference]	LPI(13:0)	1	12.4	<1
LPI(20:4) [sn2]	638.3 → 361.3	LPI 13:0 (IS) [reference]	LPI(13:0)	1	11.9	<1
methyl-CE(18:0)	684.6 → 383.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1	14.6	4.15
methyl-CE(18:1)	682.6 → 383.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1	4.3	<1
methyl-CE(18:2)	680.6 → 383.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1	2.9	<1
methyl-CE(20:4)	704.6 → 383.3	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1	4.1	<1
methyl-DE(18:1)	680.6 → 381.4	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1	9.7	<1
methyl-DE(18:2)	678.6 → 381.4	CE 18:0-d6 (IS) [reference]	CE(18:0) d6	1	6.2	<1
PA(34:1)	692.6 → 577.6	PA(15:0_18:1) d7 (IS) [reference]	PA(15:0_18:1) d7	1	11.1	<1
PA(36:2)	718.6 → 603.6	PA(15:0_18:1) d7 (IS) [reference]	PA(15:0_18:1) d7	1	13.2	5.97
PA(36:3)	716.6 → 601.6	PA(15:0_18:1) d7 (IS) [reference]	PA(15:0_18:1) d7	1	13.7	6.68
PA(36:4)	714.6 → 599.6	PA(15:0_18:1) d7 (IS) [reference]	PA(15:0_18:1) d7	1	15.5	10.42
PC(14:0_16:0)	706.5 → 184.1	PC(14:0_16:0) [reference]	PC(15:0_18:1) d7	1	1.9	<1
PC(14:0_20:4)	754.5 → 184.1	PC(14:0_20:4) [reference]	PC(15:0_18:1) d7	1	2.3	<1
PC(14:0_22:6)	778.5 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	2.7	<1
PC(15:0_20:3)	770.6 → 184.1	PC(33:2) [reference]	PC(15:0_18:1) d7	1	3.7	<1
PC(15:0_20:4)	768.6 → 184.1	PC(15:0_20:4) [reference]	PC(15:0_18:1) d7	1	2.0	<1
PC(15:0_22:6)	792.6 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	2.2	<1
PC(15-MHDA_18:1)	774.6 → 184.1	PC(16:0_18:1) [reference]	PC(15:0_18:1) d7	1	2.4	<1
PC(15-MHDA_18:2)	772.6 → 184.1	PC(16:0_18:1) [reference]	PC(15:0_18:1) d7	1	2.1	<1
PC(15-MHDA_20:4)	796.6 → 184.1	PC(16:0_18:2) [reference]	PC(15:0_18:1) d7	1	2.2	<1
PC(15-MHDA_22:6)	820.6 → 184.1	PC(18:0_22:6) [reference]	PC(15:0_18:1) d7	1	2.1	<1
PC(16:0_16:0)	734.6 → 184.1	PC(16:0_16:0) [reference]	PC(15:0_18:1) d7	1	2.0	<1
PC(16:0_18:0)	762.6 → 184.1	PC(18:0_18:1) [reference]	PC(15:0_18:1) d7	1	2.4	<1
PC(16:0_18:1)	760.6 → 184.1	PC(16:0_18:1) [reference]	PC(15:0_18:1) d7	1	3.2	<1
PC(16:0_18:2)	758.6 → 184.1	PC(16:0_18:2) [reference]	PC(15:0_18:1) d7	1	5.5	<1
PC(16:0_18:3) (a)	756.6 → 184.1	PC(32:2) [reference]	PC(15:0_18:1) d7	1	1.9	<1
PC(16:0_18:3) (b)	756.6 → 184.1	PC(32:2) [reference]	PC(15:0_18:1) d7	1	2.2	<1
PC(16:0_20:3) (a)	784.6 → 184.1	PC(33:1) [reference]	PC(15:0_18:1) d7	1	2.4	<1
PC(16:0_20:3) (b)	784.6 → 184.1	PC(33:1) [reference]	PC(15:0_18:1) d7	1	3.0	<1
PC(16:0_20:4)	782.6 → 184.1	PC(33:2) [reference]	PC(15:0_18:1) d7	1	4.0	<1
PC(16:0_20:5)	780.6 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	1.5	<1
PC(16:0_22:6)	806.6 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	2.2	<1
PC(16:1_18:2)	756.6 → 184.1	PC(33:2) [reference]	PC(15:0_18:1) d7	1	2.9	<1
PC(16:1_20:4)	780.6 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	2.9	<1
PC(16:1_22:6)	804.6 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	2.5	<1
PC(17:0_18:1)	774.6 → 184.1	PC(16:0_18:1) [reference]	PC(15:0_18:1) d7	1	2.4	<1

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
PC(17:0_18:2)	772.6 → 184.1	PC(16:0_18:1) [reference]	PC(15:0_18:1) d7	1	2.6	<1
PC(17:0_20:4)	796.6 → 184.1	PC(16:0_18:2) [reference]	PC(15:0_18:1) d7	1	2.2	<1
PC(17:0_22:6)	820.6 → 184.1	PC(18:0_22:6) [reference]	PC(15:0_18:1) d7	1	2.0	<1
PC(17:1_18:2)	770.6 → 184.1	PC(33:2) [reference]	PC(15:0_18:1) d7	1	2.3	<1
PC(18:0_18:1)	788.6 → 184.1	PC(18:0_18:1) [reference]	PC(15:0_18:1) d7	1	1.6	<1
PC(18:0_18:2)	786.6 → 184.1	PC(16:0_18:1) [reference]	PC(15:0_18:1) d7	1	3.3	<1
PC(18:0_20:3)	812.6 → 184.1	PC(18:0_18:1) [reference]	PC(15:0_18:1) d7	1	1.8	<1
PC(18:0_20:4)	810.6 → 184.1	PC(18:0_20:4) [reference]	PC(15:0_18:1) d7	1	2.0	<1
PC(18:0_22:4)	838.6 → 184.1	PC(15:0_18:1) d7 (IS) [reference]	PC(15:0_18:1) d7	1	1.6	<1
PC(18:0_22:5) (n3)/PC(20:1_20:4)	836.6 → 184.1	PC(15:0_18:1) d7 (IS) [reference]	PC(15:0_18:1) d7	1	1.8	<1
PC(18:0_22:5) (n6)	836.6 → 184.1	PC(15:0_18:1) d7 (IS) [reference]	PC(15:0_18:1) d7	1	2.1	<1
PC(18:0_22:6)	834.6 → 184.1	PC(18:0_22:6) [reference]	PC(15:0_18:1) d7	1	1.6	<1
PC(18:1_18:1)	786.6 → 184.1	PC(16:0_18:1) [reference]	PC(15:0_18:1) d7	1	3.3	<1
PC(18:1_18:2)	784.6 → 184.1	PC(33:1) [reference]	PC(15:0_18:1) d7	1	1.9	<1
PC(18:1_20:3)	810.6 → 184.1	PC(18:0_20:4) [reference]	PC(15:0_18:1) d7	1	1.6	<1
PC(18:1_22:6) (a)	832.6 → 184.1	PC(18:0_22:6) [reference]	PC(15:0_18:1) d7	1	1.6	<1
PC(18:1_22:6) (b)	832.6 → 184.1	PC(18:0_22:6) [reference]	PC(15:0_18:1) d7	1	3.9	<1
PC(18:2_18:2)	782.6 → 184.1	PC(33:2) [reference]	PC(15:0_18:1) d7	1	3.7	<1
PC(18:2_20:5)	804.6 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	2.0	<1
PC(20:0_20:4)	838.6 → 184.1	PC(15:0_18:1) d7 (IS) [reference]	PC(15:0_18:1) d7	1	3.1	<1
PC(28:0)	678.5 → 184.1	PC(28:0) [reference]	PC(15:0_18:1) d7	1	3.5	<1
PC(31:0) (a)	720.6 → 184.1	PC(16:0_18:1) [reference]	PC(15:0_18:1) d7	1	2.0	<1
PC(31:0) (b)	720.6 → 184.1	PC(16:0_18:1) [reference]	PC(15:0_18:1) d7	1	2.9	<1
PC(31:1)	718.5 → 184.1	PC(14:0_16:0) [reference]	PC(15:0_18:1) d7	1	2.4	<1
PC(32:1)	732.6 → 184.1	PC(16:0_16:0) [reference]	PC(15:0_18:1) d7	1	1.6	<1
PC(32:2)	730.5 → 184.1	PC(32:2) [reference]	PC(15:0_18:1) d7	1	1.8	<1
PC(33:0) (a)	748.6 → 184.1	PC(16:0_18:1) [reference]	PC(15:0_18:1) d7	1	1.7	<1
PC(33:0) (b)	748.6 → 184.1	PC(16:0_18:1) [reference]	PC(15:0_18:1) d7	1	2.6	<1
PC(33:1)	746.6 → 184.1	PC(33:1) [reference]	PC(15:0_18:1) d7	1	1.7	<1
PC(33:2)	744.6 → 184.1	PC(33:2) [reference]	PC(15:0_18:1) d7	1	2.8	<1
PC(34:2) [+OH]	774.6 → 184.1	PC(15:0_18:1) d7 (IS) [reference]	PC(15:0_18:1) d7	1	50.5	<1
PC(34:5)	752.5 → 184.1	PC(34:5) [reference]	PC(15:0_18:1) d7	1	3.6	<1
PC(35:5)	766.5 → 184.1	PC(15:0_20:4) [reference]	PC(15:0_18:1) d7	1	3.1	<1
PC(36:4) [+OH]	798.6 → 184.1	PC(15:0_18:1) d7 (IS) [reference]	PC(15:0_18:1) d7	1	4.2	<1
PC(36:6) (a)	778.5 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	3.0	<1
PC(38:2)	814.6 → 184.1	PC(18:0_18:1) [reference]	PC(15:0_18:1) d7	1	2.4	2.16
PC(38:4) (b)	810.6 → 184.1	PC(18:0_20:4) [reference]	PC(15:0_18:1) d7	1	1.8	<1
PC(38:5) (a)	808.6 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	2.2	<1
PC(38:5) (b)	808.6 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	1.8	<1
PC(38:6) (a)	806.6 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	1.4	<1
PC(38:6) [+OH]	822.6 → 184.1	PC(18:0_20:4) [reference]	PC(15:0_18:1) d7	1	4.3	<1
PC(38:7) (c)	804.6 → 184.1	PC(15:0_22:6) [reference]	PC(15:0_18:1) d7	1	3.3	<1
PC(39:5) (a)	822.6 → 184.1	PC(18:0_22:6) [reference]	PC(15:0_18:1) d7	1	3.1	<1
PC(39:5) (b)	822.6 → 184.1	PC(18:0_22:6) [reference]	PC(15:0_18:1) d7	1	1.9	<1
PC(40:7) (a)	832.6 → 184.1	PC(18:0_22:6) [reference]	PC(15:0_18:1) d7	1	2.7	<1
PC(40:8)	830.6 → 184.1	PC(16:0_18:2) [reference]	PC(15:0_18:1) d7	1	2.5	<1
PC(42:10)	854.5 → 184.1	PC(34:5) [reference]	PC(15:0_18:1) d7	1	3.0	<1

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
PC(O-16:0/16:0)	720.6 → 184.1	PC(O-16:0/16:0) [reference]	PC(15:0_18:1) d7	1	2.7	<1
PC(O-16:0/20:3)	770.6 → 184.1	PC(O-38:5) [reference]	PC(15:0_18:1) d7	1	2.1	<1
PC(O-16:0/20:4)	768.6 → 184.1	PC(O-38:5) [reference]	PC(15:0_18:1) d7	1	1.6	<1
PC(O-16:0/22:6)	792.6 → 184.1	PC(O-38:5) [reference]	PC(15:0_18:1) d7	1	1.6	<1
PC(O-18:0/18:1)	774.6 → 184.1	PC(O-36:0) [reference]	PC(15:0_18:1) d7	1	4.6	<1
PC(O-18:0/18:2)	772.6 → 184.1	PC(O-34:1) [reference]	PC(15:0_18:1) d7	1	2.3	<1
PC(O-18:0/20:4)	796.6 → 184.1	PC(O-38:5) [reference]	PC(15:0_18:1) d7	1	2.4	<1
PC(O-18:0/22:6)	820.6 → 184.1	PC(O-38:5) [reference]	PC(15:0_18:1) d7	1	2.6	<1
PC(O-18:1/18:1)	772.6 → 184.1	PC(O-34:1) [reference]	PC(15:0_18:1) d7	1	2.6	<1
PC(O-18:1/18:2)	770.6 → 184.1	PC(O-38:5) [reference]	PC(15:0_18:1) d7	1	1.6	<1
PC(O-32:1)	718.5 → 184.1	PC(O-16:0/16:0) [reference]	PC(15:0_18:1) d7	1	2.8	<1
PC(O-32:2)	716.6 → 184.1	PC(16:0_18:2) [reference]	PC(15:0_18:1) d7	1	2.5	1.55
PC(O-34:1)	746.6 → 184.1	PC(O-34:1) [reference]	PC(15:0_18:1) d7	1	2.0	<1
PC(O-34:2)	744.6 → 184.1	PC(O-34:1) [reference]	PC(15:0_18:1) d7	1	1.7	<1
PC(O-34:4)	740.6 → 184.1	PC(O-34:1) [reference]	PC(15:0_18:1) d7	1	2.8	2.57
PC(O-35:4)	754.5 → 184.1	PC(O-34:1) [reference]	PC(15:0_18:1) d7	1	1.9	93.31
PC(O-36:0)	776.6 → 184.1	PC(O-36:0) [reference]	PC(15:0_18:1) d7	1	3.7	<1
PC(O-36:5)	766.5 → 184.1	PC(O-38:5) [reference]	PC(15:0_18:1) d7	1	1.9	<1
PC(O-38:5)	794.6 → 184.1	PC(O-38:5) [reference]	PC(15:0_18:1) d7	1	1.9	<1
PC(O-40:5)	822.6 → 184.1	PC(O-38:5) [reference]	PC(15:0_18:1) d7	1	2.1	<1
PC(O-40:7)	818.6 → 184.1	PC(O-38:5) [reference]	PC(15:0_18:1) d7	1	2.4	<1
PC(P-15:0/20:4) (a)	752.6 → 184.1	PC(P-16:0/22:6) [reference]	PC(P-18:0/18:1) d9	1	4.5	2.73
PC(P-15:0/20:4) (b)	752.6 → 184.1	PC(P-16:0/22:6) [reference]	PC(P-18:0/18:1) d9	1	4.9	7.07
PC(P-16:0/14:0)	690.4 → 184.1	PC(P-16:0/18:1) [reference]	PC(P-18:0/18:1) d9	1	4.0	<1
PC(P-16:0/16:0)	718.5 → 184.1	PC(P-16:0/18:1) [reference]	PC(P-18:0/18:1) d9	1	3.3	<1
PC(P-16:0/16:1)	716.6 → 184.1	PC(P-16:0/18:1) [reference]	PC(P-18:0/18:1) d9	1	2.6	<1
PC(P-16:0/18:0)	746.6 → 184.1	PC(P-18:0/18:1) d9 (IS) [reference]	PC(P-18:0/18:1) d9	1	2.7	3.56
PC(P-16:0/18:1)	744.6 → 184.1	PC(P-16:0/18:1) [reference]	PC(P-18:0/18:1) d9	1	1.2	<1
PC(P-16:0/18:2)	742.5 → 184.1	PC(P-16:0/18:1) [reference]	PC(P-18:0/18:1) d9	1	3.4	<1
PC(P-16:0/18:3)	740.6 → 184.1	PC(P-16:0/20:4) [reference]	PC(P-18:0/18:1) d9	1	3.1	1.53
PC(P-16:0/20:4)	766.5 → 184.1	PC(P-16:0/20:4) [reference]	PC(P-18:0/18:1) d9	1	2.4	<1
PC(P-16:0/20:5)	764.6 → 184.1	PC(P-16:0/20:4) [reference]	PC(P-18:0/18:1) d9	1	3.3	<1
PC(P-16:0/22:6)	790.6 → 184.1	PC(P-16:0/22:6) [reference]	PC(P-18:0/18:1) d9	1	3.5	<1
PC(P-17:0/20:4) (a)	780.5 → 184.1	PC(P-16:0/20:4) [reference]	PC(P-18:0/18:1) d9	1	3.2	<1
PC(P-17:0/20:4) (b)	780.5 → 184.1	PC(P-16:0/20:4) [reference]	PC(P-18:0/18:1) d9	1	3.8	<1
PC(P-18:0/18:2)	770.6 → 184.1	PC(P-16:0/18:1) [reference]	PC(P-18:0/18:1) d9	1	1.4	<1
PC(P-18:0/20:4)	794.6 → 184.1	PC(P-16:0/20:4) [reference]	PC(P-18:0/18:1) d9	1	1.4	<1
PC(P-18:0/22:5)	820.6 → 184.1	PC(P-16:0/22:6) [reference]	PC(P-18:0/18:1) d9	1	1.8	<1
PC(P-18:0/22:6)	818.6 → 184.1	PC(P-16:0/22:6) [reference]	PC(P-18:0/18:1) d9	1	2.9	<1
PC(P-18:1/18:1)	770.6 → 184.1	PC(P-16:0/18:1) [reference]	PC(P-18:0/18:1) d9	1	2.1	<1
PC(P-18:1/22:6)	816.6 → 184.1	PC(P-16:0/22:6) [reference]	PC(P-18:0/18:1) d9	1	5.8	1.18
PC(P-20:0/20:4)	822.6 → 184.1	PC(P-18:0/18:1) d9 (IS) [reference]	PC(P-18:0/18:1) d9	1	2.6	1.89
PC(P-35:2) (a)	756.6 → 184.1	PC(P-16:0/18:1) [reference]	PC(P-18:0/18:1) d9	1	2.9	<1
PC(P-35:2) (b)	756.6 → 184.1	PC(P-16:0/18:1) [reference]	PC(P-18:0/18:1) d9	1	3.0	<1
PC(P-36:3)	768.5 → 184.1	PC(P-16:0/20:4) [reference]	PC(P-18:0/18:1) d9	1	2.1	<1
PC(P-38:5) (a)	792.6 → 184.1	PC(P-16:0/22:6) [reference]	PC(P-18:0/18:1) d9	1	3.1	<1
PC(P-38:5) (b)	792.6 → 184.1	PC(P-16:0/22:6) [reference]	PC(P-18:0/18:1) d9	1	2.2	<1

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
PE(15-MHDA_18:1)	732.6 → 591.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	12.8	17.5
PE(15-MHDA_18:2)	730.5 → 589.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	9.1	<1
PE(15-MHDA_20:4)	754.6 → 613.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	11.1	<1
PE(15-MHDA_22:6)	778.5 → 637.5	PE(16:0_22:6) [reference]	PE(15:0_18:1) d7	1	9.4	<1
PE(16:0_16:0)	692.5 → 551.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	17.5	<1
PE(16:0_16:1)	690.5 → 549.5	PE(16:0_16:1) [reference]	PE(15:0_18:1) d7	1	8.9	<1
PE(16:0_18:1)	718.5 → 577.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	2.5	<1
PE(16:0_18:2)	716.5 → 575.5	PE(16:0_18:2) [reference]	PE(15:0_18:1) d7	1	2.5	<1
PE(16:0_18:3) (a)	714.5 → 573.5	PE(16:0_18:2) [reference]	PE(15:0_18:1) d7	1	7.0	1.14
PE(16:0_18:3) (b)	714.5 → 573.5	PE(16:0_18:2) [reference]	PE(15:0_18:1) d7	1	6.9	<1
PE(16:0_20:3)	742.5 → 601.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	2.7	<1
PE(16:0_20:4)	740.5 → 599.5	PE(18:0_20:4) [reference]	PE(15:0_18:1) d7	1	1.5	<1
PE(16:0_20:5)	738.5 → 597.5	PE(16:0_16:1) [reference]	PE(15:0_18:1) d7	1	4.3	<1
PE(16:0_22:6)	764.5 → 623.5	PE(16:0_22:6) [reference]	PE(15:0_18:1) d7	1	5.8	<1
PE(16:1_18:2)	714.5 → 573.5	PE(16:0_18:2) [reference]	PE(15:0_18:1) d7	1	10.5	<1
PE(16:1_20:4)	738.5 → 597.5	PE(16:0_16:1) [reference]	PE(15:0_18:1) d7	1	7.3	<1
PE(17:0_18:1)	732.6 → 591.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	11.0	<1
PE(17:0_18:2)	730.5 → 589.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	8.0	1.13
PE(17:0_20:4)	754.6 → 613.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	5.0	<1
PE(17:0_22:6)	778.5 → 637.5	PE(16:0_22:6) [reference]	PE(15:0_18:1) d7	1	9.1	2.42
PE(18:0_18:1)	746.6 → 605.6	PE(18:0_18:1) [reference]	PE(15:0_18:1) d7	1	2.0	<1
PE(18:0_18:2)	744.6 → 603.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	1.5	<1
PE(18:0_20:3) (a)	770.6 → 629.6	PE(18:0_18:1) [reference]	PE(15:0_18:1) d7	1	2.3	<1
PE(18:0_20:3) (b)	770.6 → 629.6	PE(18:0_18:1) [reference]	PE(15:0_18:1) d7	1	4.3	<1
PE(18:0_20:4)	768.6 → 627.5	PE(18:0_20:4) [reference]	PE(15:0_18:1) d7	1	1.4	<1
PE(18:0_22:4)	796.6 → 655.6	PE(18:0_18:1) [reference]	PE(15:0_18:1) d7	1	4.5	<1
PE(18:0_22:5) (n3)	794.6 → 653.6	PE(18:0_18:1) [reference]	PE(15:0_18:1) d7	1	2.5	<1
PE(18:0_22:5) (n6)	794.6 → 653.6	PE(18:0_18:1) [reference]	PE(15:0_18:1) d7	1	3.8	<1
PE(18:0_22:6)	792.6 → 651.5	PE(18:0_22:6) [reference]	PE(15:0_18:1) d7	1	2.9	<1
PE(18:1_18:1)	744.6 → 603.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	5.8	<1
PE(18:1_18:2)	742.5 → 601.5	PE(16:0_18:1) [reference]	PE(15:0_18:1) d7	1	3.0	<1
PE(18:1_22:6) (a)	790.5 → 649.5	PE(18:0_22:6) [reference]	PE(15:0_18:1) d7	1	4.2	1.05
PE(18:1_22:6) (b)	790.5 → 649.5	PE(18:0_22:6) [reference]	PE(15:0_18:1) d7	1	7.9	1.59
PE(20:0_20:4)	796.6 → 655.6	PE(18:0_18:1) [reference]	PE(15:0_18:1) d7	1	7.8	<1
PE(38:5) (a)	766.5 → 625.5	PE(18:0_20:4) [reference]	PE(15:0_18:1) d7	1	3.2	<1
PE(38:5) (b)	766.5 → 625.5	PE(18:0_20:4) [reference]	PE(15:0_18:1) d7	1	2.7	<1
PE(O-16:0/18:2)	702.5 → 561.5	PE(O-34:1) [reference]	PE(15:0_18:1) d7	1	4.6	<1
PE(O-16:0/20:3)	728.6 → 587.5	PE(18:0_22:6) [reference]	PE(15:0_18:1) d7	1	6.9	<1
PE(O-16:0/20:4)	726.5 → 585.5	PE(P-18:0/20:4) [reference]	PE(15:0_18:1) d7	1	2.5	<1
PE(O-16:0/22:4)	754.6 → 613.6	PE(P-18:0/20:4) [reference]	PE(15:0_18:1) d7	1	5.1	8.97
PE(O-16:0/22:6)	750.6 → 609.5	PE(P-18:0/20:4) [reference]	PE(15:0_18:1) d7	1	2.5	<1
PE(O-18:0/20:4)	754.6 → 613.6	PE(P-18:0/20:4) [reference]	PE(15:0_18:1) d7	1	2.7	<1
PE(O-18:0/22:5)	780.6 → 639.6	PE(P-18:0/20:4) [reference]	PE(15:0_18:1) d7	1	5.1	<1
PE(O-18:0/22:6)	778.5 → 637.5	PE(P-18:0/20:4) [reference]	PE(15:0_18:1) d7	1	3.1	<1
PE(O-18:1/18:2)	728.6 → 587.5	PE(18:0_22:6) [reference]	PE(15:0_18:1) d7	1	4.5	<1
PE(O-18:1/22:6)	776.6 → 635.5	PE(P-18:0/20:4) [reference]	PE(15:0_18:1) d7	1	2.9	<1
PE(O-34:1)	704.6 → 563.5	PE(O-34:1) [reference]	PE(15:0_18:1) d7	1	3.1	<1

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
PE(O-36:5)	724.5 → 583.5	PE(P-18:0/20:4) [reference]	PE(15:0_18:1) d7	1	6.5	1.71
PE(O-38:5) (a)	752.6 → 611.5	PE(P-18:0/20:4) [reference]	PE(15:0_18:1) d7	1	3.0	<1
PE(O-38:5) (b)	752.6 → 611.5	PE(P-18:0/20:4) [reference]	PE(15:0_18:1) d7	1	4.7	1.69
PE(P-15:0/22:6)	734.5 → 385.3	PE(P-16:0/18:2) [reference]	PE(P-18:0/18:1) d9	1	12.0	<1
PE(P-16:0/18:1)	702.5 → 339.3	PE(P-16:0/18:2) [reference]	PE(P-18:0/18:1) d9	1	2.8	<1
PE(P-16:0/18:2)	700.5 → 337.3	PE(P-16:0/18:2) [reference]	PE(P-18:0/18:1) d9	1	2.1	<1
PE(P-16:0/20:3) (a)	726.5 → 363.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	4.9	<1
PE(P-16:0/20:3) (b)	726.5 → 363.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	8.0	1.32
PE(P-16:0/20:4)	724.5 → 361.3	PE(P-16:0/18:2) [reference]	PE(P-18:0/18:1) d9	1	3.0	<1
PE(P-16:0/20:5)	722.5 → 359.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	5.2	<1
PE(P-16:0/22:4)	752.6 → 389.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	5.2	<1
PE(P-16:0/22:5) (n3)	750.5 → 387.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	2.6	<1
PE(P-16:0/22:5) (n6)	750.5 → 387.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	6.2	<1
PE(P-16:0/22:6)	748.5 → 385.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	2.5	<1
PE(P-17:0/20:4) (a)	738.6 → 361.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	4.7	<1
PE(P-17:0/20:4) (b)	738.6 → 361.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	5.2	<1
PE(P-17:0/22:6) (a)	762.6 → 385.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	5.9	<1
PE(P-17:0/22:6) (b)	762.6 → 385.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	5.9	<1
PE(P-18:0/18:1)	730.6 → 339.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	2.9	<1
PE(P-18:0/18:2)	728.6 → 337.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	2.1	<1
PE(P-18:0/18:3)	726.5 → 335.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	12.7	<1
PE(P-18:0/20:3) (a)	754.5 → 363.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	3.6	<1
PE(P-18:0/20:3) (b)	754.5 → 363.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	6.8	<1
PE(P-18:0/20:4)	752.6 → 361.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	1.4	<1
PE(P-18:0/20:5)	750.5 → 359.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	2.8	<1
PE(P-18:0/22:4)	780.6 → 389.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	4.9	<1
PE(P-18:0/22:5) (n3)	778.5 → 387.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	3.2	<1
PE(P-18:0/22:5) (n6)	778.5 → 387.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	6.7	1.39
PE(P-18:0/22:6)	776.6 → 385.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	2.0	<1
PE(P-18:1/18:1) (a)	728.6 → 339.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	3.0	<1
PE(P-18:1/18:1) (b)	728.6 → 339.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	6.2	<1
PE(P-18:1/18:2) (a)	726.5 → 337.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	3.0	<1
PE(P-18:1/18:2) (b)	726.5 → 337.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	5.5	<1
PE(P-18:1/20:3) (a)	752.5 → 363.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	6.1	<1
PE(P-18:1/20:3) (b)	752.5 → 363.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	9.5	<1
PE(P-18:1/20:4) (a)	750.5 → 361.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	2.4	<1
PE(P-18:1/20:4) (b)	750.5 → 361.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	5.0	<1
PE(P-18:1/20:5) (a)	748.5 → 359.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	4.1	<1
PE(P-18:1/20:5) (b)	748.5 → 359.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	7.7	<1
PE(P-18:1/22:4)	778.5 → 389.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	9.4	<1
PE(P-18:1/22:5) (a)	776.6 → 387.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	5.4	2.73
PE(P-18:1/22:5) (b)	776.6 → 387.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	8.9	<1
PE(P-18:1/22:6) (a)	774.5 → 385.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	1.9	<1
PE(P-18:1/22:6) (b)	774.5 → 385.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	4.1	<1
PE(P-19:0/20:4) (a)	766.6 → 361.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	12.3	1.51
PE(P-19:0/20:4) (b)	766.6 → 361.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	9.2	<1
PE(P-20:0/18:1)	758.6 → 339.3	PE(P-18:0/18:1) d9 (IS) [reference]	PE(P-18:0/18:1) d9	1	6.7	3.32



化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
PE(P-20:0/18:2)	756.6 → 337.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	4.7	<1
PE(P-20:0/20:4)	780.6 → 361.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	3.3	<1
PE(P-20:0/22:6)	804.6 → 385.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	4.0	<1
PE(P-20:1/20:4)	778.5 → 361.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	3.8	<1
PE(P-20:1/22:6)	802.6 → 385.3	PE(P-18:0/20:4) [reference]	PE(P-18:0/18:1) d9	1	5.5	<1
PG(34:1)	766.6 → 577.5	PG(15:0_18:1) d7 (IS) [reference]	PG(15:0_18:1) d7	1	12.2	<1
PG(36:1)	794.6 → 605.6	PG(15:0_18:1) d7 (IS) [reference]	PG(15:0_18:1) d7	1	10.9	<1
PG(36:2)	792.6 → 603.5	PG(15:0_18:1) d7 (IS) [reference]	PG(15:0_18:1) d7	1	12.2	<1
PI(38:5) (b)	902.6 → 625.6	PI(38:6) [reference]	PI(15:0_18:1) d7	1	4.0	<1
PI(15-MHDA_18:1)/PI(17:0_18:1)	868.6 → 591.6	PI(34:1) [reference]	PI(15:0_18:1) d7	1	7.3	<1
PI(15-MHDA_18:2)/PI(17:0_18:2)	866.6 → 589.6	PI(15-MHDA_18:2)/PI(17:0_18:2) [reference]	PI(15:0_18:1) d7	1	8.0	<1
PI(37:4)	890.6 → 613.6	PI(16:0_20:4) [reference]	PI(15:0_18:1) d7	1	6.4	<1
PI(16:0/16:0)	828.6 → 551.6	PI(34:1) [reference]	PI(15:0_18:1) d7	1	5.7	<1
PI(16:0_16:1)	826.5 → 549.5	PI(16:0_16:1) [reference]	PI(15:0_18:1) d7	1	3.8	<1
PI(16:0_20:3) (a)	878.6 → 601.6	PI(16:0_20:4) [reference]	PI(15:0_18:1) d7	1	4.7	<1
PI(16:0_20:3) (b)	878.6 → 601.6	PI(16:0_20:4) [reference]	PI(15:0_18:1) d7	1	5.3	<1
PI(16:0_20:4)	876.6 → 599.6	PI(16:0_20:4) [reference]	PI(15:0_18:1) d7	1	2.2	<1
PI(18:0_18:1)	882.6 → 605.6	PI(18:0_18:1) [reference]	PI(15:0_18:1) d7	1	2.5	<1
PI(18:0_20:2)	908.6 → 631.6	PI(18:0_20:4) [reference]	PI(15:0_18:1) d7	1	5.7	<1
PI(18:0_20:3) (a)	906.6 → 629.6	PI(18:0_20:4) [reference]	PI(15:0_18:1) d7	1	9.5	<1
PI(18:0_20:3) (b)	906.6 → 629.6	PI(18:0_20:4) [reference]	PI(15:0_18:1) d7	1	3.2	<1
PI(18:0_20:4)	904.6 → 627.6	PI(18:0_20:4) [reference]	PI(15:0_18:1) d7	1	3.7	<1
PI(18:0_22:4)	932.6 → 655.6	PI(18:0_22:6) [reference]	PI(15:0_18:1) d7	1	6.2	<1
PI(18:0_22:5) (n3)	930.6 → 653.6	PI(18:0_22:6) [reference]	PI(15:0_18:1) d7	1	13.4	<1
PI(18:0_22:5) (n6)	930.6 → 653.6	PI(18:0_22:6) [reference]	PI(15:0_18:1) d7	1	9.0	<1
PI(18:0_22:6)	928.6 → 651.6	PI(18:0_22:6) [reference]	PI(15:0_18:1) d7	1	3.3	<1
PI(18:1_18:2)	878.6 → 601.6	PI(16:0_20:4) [reference]	PI(15:0_18:1) d7	1	3.8	<1
PI(20:0_20:4)	932.6 → 655.6	PI(18:0_22:6) [reference]	PI(15:0_18:1) d7	1	9.2	<1
PI(34:0)	856.6 → 579.6	PI(18:0_18:1) [reference]	PI(15:0_18:1) d7	1	7.8	<1
PI(34:1)	854.6 → 577.6	PI(34:1) [reference]	PI(15:0_18:1) d7	1	2.3	<1
PI(36:2)	880.6 → 603.6	PI(16:0_20:4) [reference]	PI(15:0_18:1) d7	1	10.9	<1
PI(37:6)	886.6 → 609.6	PI(16:0_20:4) [reference]	PI(15:0_18:1) d7	1	11.7	<1
PI(38:5) (a)	902.6 → 625.6	PI(38:6) [reference]	PI(15:0_18:1) d7	1	3.7	<1
PI(38:6)	900.6 → 623.6	PI(38:6) [reference]	PI(15:0_18:1) d7	1	3.6	<1
PIP1(38:4)	984.7 → 627.7	PI(18:0_22:6) [reference]	PI(15:0_18:1) d7	1	7.0	1.52
PS(36:1)	790.6 → 605.6	PS(15:0_18:1) d7 (IS) [reference]	PS(15:0_18:1) d7	1	4.7	<1
PS(36:2)	788.5 → 603.5	PS(15:0_18:1) d7 (IS) [reference]	PS(15:0_18:1) d7	1	4.9	6.95
PS(38:4)	812.5 → 627.5	PS(15:0_18:1) d7 (IS) [reference]	PS(15:0_18:1) d7	1	11.9	<1
PS(40:5)	838.6 → 653.6	PS(15:0_18:1) d7 (IS) [reference]	PS(15:0_18:1) d7	1	6.8	16.53
S1P(d16:1)	352.2 → 236.3	S1P(d18:1) d7 (IS) [reference]	S1P(18:1) d7	1	6.9	4.6
S1P(d18:0)	382.2 → 284.3	S1P(d18:1) d7 (IS) [reference]	S1P(18:1) d7	1	14.7	<1
S1P(d18:1)	380.2 → 264.3	S1P(d18:1) d7 (IS) [reference]	S1P(18:1) d7	1	5.6	<1
S1P(d18:2)	378.2 → 262.3	S1P(d18:1) d7 (IS) [reference]	S1P(18:1) d7	1	5.1	<1
SHexCer(d18:1/16:0(OH))	796.8 → 264.3	SHexCer(d18:1/16:0) [reference]	SHexCer(d18:1/12:0)	1	6.2	<1
SHexCer(d18:1/16:0)	780.8 → 264.3	SHexCer(d18:1/16:0) [reference]	SHexCer(d18:1/12:0)	1	8.1	<1
SHexCer(d18:1/24:0(OH))	908.8 → 264.3	SHexCer(d18:1/24:1(OH)) [reference]	SHexCer(d18:1/12:0)	1	10.9	<1
SHexCer(d18:1/24:1(OH))	906.8 → 264.3	SHexCer(d18:1/24:1(OH)) [reference]	SHexCer(d18:1/12:0)	1	12.8	<1

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
SHexCer(d18:1/24:1)	890.8 → 264.3	SHexCer(d18:1/24:1(OH)) [reference]	SHexCer(d18:1/12:0)	1	16.3	<1
SM(34:3)	699.5 → 184.1	SM(d18:2/14:0) [reference]	SM(d18:1/15:0) d9	1	4.4	<1
SM(35:2) (b)	715.6 → 184.1	SM(d17:1/16:0) [reference]	SM(d18:1/15:0) d9	1	5.0	<1
SM(37:1)	745.6 → 184.1	SM(d18:1/18:0)/SM(d16:1/20:0) [reference]	SM(d18:1/15:0) d9	1	1.8	<1
SM(37:2)	743.5 → 184.1	SM(d18:2/18:0) [reference]	SM(d18:1/15:0) d9	1	1.6	<1
SM(38:3) (a)	755.6 → 184.1	SM(d18:2/20:0) [reference]	SM(d18:1/15:0) d9	1	3.0	<1
SM(38:3) (b)	755.6 → 184.1	SM(d18:2/20:0) [reference]	SM(d18:1/15:0) d9	1	2.8	66.33
SM(40:3) (a)	783.6 → 184.1	SM(d18:2/20:0) [reference]	SM(d18:1/15:0) d9	1	2.2	<1
SM(40:3) (b)	783.6 → 184.1	SM(d18:2/20:0) [reference]	SM(d18:1/15:0) d9	1	2.0	<1
SM(40:4)	781.5 → 184.1	SM(d18:2/20:0) [reference]	SM(d18:1/15:0) d9	1	4.5	<1
SM(41:0)	803.7 → 184.1	SM(d18:1/24:0) [reference]	SM(d18:1/15:0) d9	1	3.3	<1
SM(41:1) (a)	801.7 → 184.1	SM(d16:1/23:0)/SM(d17:1/22:0) [reference]	SM(d18:1/15:0) d9	1	5.3	<1
SM(43:1)	829.7 → 184.1	SM(43:1) [reference]	SM(d18:1/15:0) d9	1	2.5	<1
SM(43:2) (b)	827.7 → 184.1	SM(d19:1/24:1) [reference]	SM(d18:1/15:0) d9	1	2.5	<1
SM(43:2) (c)	827.7 → 184.1	SM(d19:1/24:1) [reference]	SM(d18:1/15:0) d9	1	2.3	<1
SM(44:1)	843.6 → 184.1	SM(44:2) [reference]	SM(d18:1/15:0) d9	1	4.5	<1
SM(44:2)	841.6 → 184.1	SM(44:2) [reference]	SM(d18:1/15:0) d9	1	2.8	<1
SM(44:3)	839.6 → 184.1	SM(44:2) [reference]	SM(d18:1/15:0) d9	1	4.2	<1
SM(d16:1/19:0)	717.6 → 184.1	SM(d18:2/18:0) [reference]	SM(d18:1/15:0) d9	1	4.0	<1
SM(d16:1/23:0)/SM(d17:1/22:0)	773.7 → 184.1	SM(d16:1/23:0)/SM(d17:1/22:0) [reference]	SM(d18:1/15:0) d9	1	2.9	<1
SM(d16:1/24:1)	785.7 → 184.1	SM(d18:1/22:0)/SM(d16:1/24:0) [reference]	SM(d18:1/15:0) d9	1	2.0	<1
SM(d17:1/14:0)	661.5 → 184.1	SM(d18:1/15:0) d9 (IS) [reference]	SM(d18:1/15:0) d9	1	2.8	<1
SM(d17:1/16:0)	689.6 → 184.1	SM(d17:1/16:0) [reference]	SM(d18:1/15:0) d9	1	1.4	<1
SM(d17:1/24:1)	799.7 → 184.1	SM(d16:1/23:0)/SM(d17:1/22:0) [reference]	SM(d18:1/15:0) d9	1	2.9	<1
SM(d18:0/14:0)	677.6 → 184.1	SM(d18:1/15:0) d9 (IS) [reference]	SM(d18:1/15:0) d9	1	4.4	<1
SM(d18:0/16:0)	705.6 → 184.1	SM(d18:1/16:0) [reference]	SM(d18:1/15:0) d9	1	4.2	<1
SM(d18:0/22:0)	789.7 → 184.1	SM(d18:1/22:0)/SM(d16:1/24:0) [reference]	SM(d18:1/15:0) d9	1	6.6	<1
SM(d18:1/14:0)/SM(d16:1/16:0)	675.5 → 184.1	SM(d18:1/15:0) d9 (IS) [reference]	SM(d18:1/15:0) d9	1	3.6	<1
SM(d18:1/16:0)	703.6 → 184.1	SM(d18:1/16:0) [reference]	SM(d18:1/15:0) d9	1	2.8	<1
SM(d18:1/17:0)/SM(d17:1/18:0)	717.6 → 184.1	SM(d18:2/18:0) [reference]	SM(d18:1/15:0) d9	1	3.0	<1
SM(d18:1/18:0)/SM(d16:1/20:0)	731.6 → 184.1	SM(d18:1/18:0)/SM(d16:1/20:0) [reference]	SM(d18:1/15:0) d9	1	2.1	<1
SM(d18:1/20:0)/SM(d16:1/22:0)	759.6 → 184.1	SM(d16:1/23:0)/SM(d17:1/22:0) [reference]	SM(d18:1/15:0) d9	1	1.7	<1
SM(d18:1/22:0)/SM(d16:1/24:0)	787.7 → 184.1	SM(d18:1/22:0)/SM(d16:1/24:0) [reference]	SM(d18:1/15:0) d9	1	1.6	<1
SM(d18:1/23:0)/SM(d17:1/24:0)	801.7 → 184.1	SM(d16:1/23:0)/SM(d17:1/22:0) [reference]	SM(d18:1/15:0) d9	1	1.3	<1
SM(d18:1/24:0)	815.7 → 184.1	SM(d18:1/24:0) [reference]	SM(d18:1/15:0) d9	1	4.0	<1
SM(d18:1/24:1)	813.7 → 184.1	SM(d18:1/24:1) [reference]	SM(d18:1/15:0) d9	1	1.7	<1
SM(d18:2/14:0)	673.5 → 184.1	SM(d18:2/14:0) [reference]	SM(d18:1/15:0) d9	1	5.6	<1
SM(d18:2/16:0)	701.6 → 184.1	SM(d18:2/16:0) [reference]	SM(d18:1/15:0) d9	1	2.8	<1
SM(d18:2/17:0)	715.6 → 184.1	SM(d17:1/16:0) [reference]	SM(d18:1/15:0) d9	1	2.9	<1
SM(d18:2/18:0)	729.6 → 184.1	SM(d18:2/18:0) [reference]	SM(d18:1/15:0) d9	1	1.9	<1
SM(d18:2/18:1)	727.6 → 184.1	SM(d18:2/16:0) [reference]	SM(d18:1/15:0) d9	1	2.2	<1
SM(d18:2/20:0)	757.6 → 184.1	SM(d18:2/20:0) [reference]	SM(d18:1/15:0) d9	1	2.0	<1
SM(d18:2/22:0)	785.7 → 184.1	SM(d18:1/22:0)/SM(d16:1/24:0) [reference]	SM(d18:1/15:0) d9	1	1.7	<1
SM(d18:2/23:0)	799.7 → 184.1	SM(d16:1/23:0)/SM(d17:1/22:0) [reference]	SM(d18:1/15:0) d9	1	2.7	<1
SM(d18:2/24:0)	813.7 → 184.1	SM(d18:1/24:1) [reference]	SM(d18:1/15:0) d9	1	4.1	<1
SM(d19:1/24:1)	827.7 → 184.1	SM(d19:1/24:1) [reference]	SM(d18:1/15:0) d9	1	4.2	<1
Sph(d18:1)	300.3 → 282.3	Sph(d17:1) [reference]	Sph(d17:1)	1	6.8	58.42

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
Sph(d18:2)	298.3 → 280.3	Sph(d17:1) [reference]	Sph(d17:1)	1	25.7	76.52
TG(48:0) [NL-16:0]	824.8 → 551.5	TG(48:0) [NL-16:0] [reference]	TG(15:0_18:1-d7_15:0)	1	4.0	1.25
TG(48:0) [NL-18:0]	824.8 → 523.5	TG(48:0) [NL-16:0] [reference]	TG(15:0_18:1-d7_15:0)	1	3.9	2.08
TG(48:0) [SIM]	824.8 → 824.8	TG(48:0) [NL-16:0] [reference]	TG(15:0_18:1-d7_15:0)	1	2.5	2.86
TG(48:1) [NL-16:1]	822.8 → 551.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.9	2.66
TG(48:1) [NL-18:1]	822.8 → 523.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.0	<1
TG(48:1) [SIM]	822.8 → 822.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	2.5	1.49
TG(48:2) [NL-14:0]	820.8 → 575.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.0	<1
TG(48:2) [NL-14:1]	820.8 → 577.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.7	1.14
TG(48:2) [NL-16:1]	820.8 → 549.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	2.7	4.14
TG(48:2) [NL-18:2]	820.8 → 523.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	2.7	<1
TG(48:2) [SIM]	820.8 → 820.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.3	<1
TG(48:3) [NL-14:0]	818.8 → 573.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.7	<1
TG(48:3) [NL-16:1]	818.8 → 547.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.3	5.58
TG(48:3) [NL-18:3]	818.8 → 523.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	4.6	<1
TG(48:3) [SIM]	818.8 → 818.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.8	4.47
TG(49:1) [NL-16:1]	836.8 → 565.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	4.9	14.31
TG(49:1) [NL-17:1]	836.8 → 551.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	2.8	<1
TG(49:1) [SIM]	836.8 → 836.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	2.3	13.44
TG(50:0) [NL-18:0]	852.8 → 551.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	4.8	<1
TG(50:0) [SIM]	852.8 → 852.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	2.3	1.84
TG(50:1) [NL-14:0]	850.8 → 605.5	TG(50:1) [NL-16:0] [reference]	TG(15:0_18:1-d7_15:0)	1	5.4	<1
TG(50:1) [NL-16:0]	850.8 → 577.5	TG(50:1) [NL-16:0] [reference]	TG(15:0_18:1-d7_15:0)	1	3.2	<1
TG(50:1) [NL-18:1]	850.8 → 551.5	TG(50:1) [NL-16:0] [reference]	TG(15:0_18:1-d7_15:0)	1	3.5	<1
TG(50:1) [SIM]	850.8 → 850.8	TG(50:1) [NL-16:0] [reference]	TG(15:0_18:1-d7_15:0)	1	3.5	<1
TG(50:2) [NL-14:0]	848.8 → 603.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	4.0	<1
TG(50:2) [NL-16:1]	848.8 → 577.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.6	<1
TG(50:2) [NL-18:1]	848.8 → 549.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.0	<1
TG(50:2) [NL-18:2]	848.8 → 551.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	4.0	<1
TG(50:2) [SIM]	848.8 → 848.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.6	<1
TG(50:3) [NL-14:0]	846.8 → 601.5	TG(50:3) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	2.9	<1
TG(50:3) [NL-14:1]	846.8 → 603.5	TG(50:3) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	4.0	<1
TG(50:3) [NL-16:1]	846.8 → 575.5	TG(50:3) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	2.4	<1
TG(50:3) [NL-18:2]	846.8 → 549.5	TG(50:3) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	2.4	<1
TG(50:3) [NL-18:3]	846.8 → 551.5	TG(50:3) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	3.1	<1
TG(50:3) [SIM]	846.8 → 846.8	TG(50:3) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	2.5	<1
TG(50:4) [NL-14:0]	844.8 → 599.5	TG(50:4) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.3	<1
TG(50:4) [NL-18:3]	844.8 → 549.5	TG(50:4) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.2	<1
TG(50:4) [NL-20:4]	844.8 → 523.5	TG(50:4) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	4.2	<1
TG(50:4) [SIM]	844.8 → 844.8	TG(50:4) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.4	<1
TG(51:0) [NL-16:0]	866.7 → 593.4	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.2	<1
TG(51:0) [SIM]	866.7 → 866.7	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	9.6	<1
TG(51:1) [NL-17:0]	864.8 → 577.5	TG(51:1) [NL-17:0] [reference]	TG(15:0_18:1-d7_15:0)	1	5.6	<1
TG(51:1) [SIM]	864.8 → 864.8	TG(51:2) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.8	<1
TG(51:2) [NL-15:0]	862.8 → 603.5	TG(51:2) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	5.6	<1
TG(51:2) [NL-17:0]	862.8 → 575.5	TG(51:2) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.9	<1
TG(51:2) [NL-17:1]	862.8 → 577.5	TG(51:2) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.3	<1

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
TG(51:2) [SIM]	862.8 → 862.8	TG(51:2) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.4	3.13
TG(52:1) [NL-18:0]	878.8 → 577.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	4.2	<1
TG(52:1) [NL-18:1]	878.8 → 579.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.3	<1
TG(52:1) [SIM]	878.8 → 878.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.5	<1
TG(52:2) [NL-16:0]	876.8 → 603.5	TG(52:2) [NL-16:0] [reference]	TG(15:0_18:1-d7_15:0)	1	4.8	<1
TG(52:2) [NL-18:2]	876.8 → 579.5	TG(52:2) [NL-16:0] [reference]	TG(15:0_18:1-d7_15:0)	1	2.8	<1
TG(52:2) [SIM]	876.8 → 876.8	TG(52:2) [NL-16:0] [reference]	TG(15:0_18:1-d7_15:0)	1	8.1	<1
TG(52:3) [NL-16:1]	874.8 → 603.5	TG(52:3) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	6.1	<1
TG(52:3) [NL-18:2]	874.8 → 577.5	TG(52:3) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	3.3	<1
TG(52:3) [SIM]	874.8 → 874.8	TG(52:3) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	8.1	<1
TG(52:4) [NL-16:1]	872.8 → 601.5	TG(52:4) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.2	<1
TG(52:4) [NL-18:2]	872.8 → 575.5	TG(52:4) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.3	<1
TG(52:4) [NL-18:3]	872.8 → 577.5	TG(52:4) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.6	<1
TG(52:4) [SIM]	872.8 → 872.8	TG(52:4) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.6	<1
TG(52:5) [NL-18:3]	870.8 → 575.5	TG(52:5) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.9	<1
TG(52:5) [NL-20:4]	870.8 → 549.5	TG(52:5) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.3	<1
TG(52:5) [NL-20:5]	870.8 → 551.5	TG(52:5) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	4.0	<1
TG(52:5) [SIM]	870.8 → 870.8	TG(52:5) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.7	<1
TG(53:2) [NL-17:1]	890.8 → 605.5	TG(53:2) [NL-18:1] [reference]	TG(15:0_18:1-d7_15:0)	1	5.5	<1
TG(53:2) [NL-18:1]	890.8 → 591.5	TG(53:2) [NL-18:1] [reference]	TG(15:0_18:1-d7_15:0)	1	2.3	<1
TG(53:2) [SIM]	890.8 → 890.8	TG(53:2) [NL-18:1] [reference]	TG(15:0_18:1-d7_15:0)	1	2.3	<1
TG(54:0) [NL-18:0]	908.8 → 607.5	TG(53:2) [NL-18:1] [reference]	TG(15:0_18:1-d7_15:0)	1	2.4	8.23
TG(54:0) [SIM]	908.8 → 908.8	TG(53:2) [NL-18:1] [reference]	TG(15:0_18:1-d7_15:0)	1	2.0	4.98
TG(54:1) [NL-18:1]	906.8 → 607.5	TG(53:2) [NL-18:1] [reference]	TG(15:0_18:1-d7_15:0)	1	2.4	<1
TG(54:1) [SIM]	906.8 → 906.8	TG(53:2) [NL-18:1] [reference]	TG(15:0_18:1-d7_15:0)	1	2.6	<1
TG(54:2) [NL-18:0]	904.8 → 603.5	TG(54:2) [NL-20:1] [reference]	TG(15:0_18:1-d7_15:0)	1	4.1	<1
TG(54:2) [NL-20:1]	904.8 → 577.5	TG(54:2) [NL-20:1] [reference]	TG(15:0_18:1-d7_15:0)	1	4.4	<1
TG(54:2) [SIM]	904.8 → 904.8	TG(54:2) [NL-20:1] [reference]	TG(15:0_18:1-d7_15:0)	1	3.7	<1
TG(54:3) [NL-18:1]	902.8 → 603.5	TG(54:3) [NL-18:1] [reference]	TG(15:0_18:1-d7_15:0)	1	2.8	<1
TG(54:3) [NL-18:2]	902.8 → 605.5	TG(54:3) [NL-18:1] [reference]	TG(15:0_18:1-d7_15:0)	1	5.4	<1
TG(54:3) [SIM]	902.8 → 902.8	TG(54:3) [NL-18:1] [reference]	TG(15:0_18:1-d7_15:0)	1	2.7	<1
TG(54:4) [NL-18:2]	900.8 → 603.5	TG(54:4) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	4.3	<1
TG(54:4) [NL-20:3]	900.8 → 577.5	TG(54:4) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	3.9	<1
TG(54:4) [SIM]	900.8 → 900.8	TG(54:4) [NL-18:2] [reference]	TG(15:0_18:1-d7_15:0)	1	3.0	<1
TG(54:5) [NL-18:3]	898.8 → 603.5	TG(54:5) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.6	<1
TG(54:5) [NL-20:4]	898.8 → 577.5	TG(54:5) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.3	<1
TG(54:5) [SIM]	898.8 → 898.8	TG(54:5) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.4	<1
TG(54:6) [NL-18:3]	896.8 → 601.5	TG(54:6) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.0	<1
TG(54:6) [NL-20:4]	896.8 → 575.5	TG(54:6) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.0	<1
TG(54:6) [NL-20:5]	896.8 → 577.5	TG(54:6) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.8	<1
TG(54:6) [NL-22:6]	896.8 → 551.5	TG(54:6) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.1	<1
TG(54:6) [SIM]	896.8 → 896.8	TG(54:6) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.2	<1
TG(54:7) [NL-20:5]	894.8 → 575.5	TG(54:7) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	3.7	<1
TG(54:7) [NL-22:6]	894.8 → 549.5	TG(54:7) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	3.5	<1
TG(54:7) [SIM]	894.8 → 894.8	TG(54:7) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	2.1	<1
TG(56:6) [NL-20:4]	924.8 → 603.5	TG(56:6) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	6.3	<1
TG(56:6) [NL-22:5]	924.8 → 577.5	TG(56:6) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.2	<1

化合物名称	离子对	时间参比化合物	内标化合物	响应因子	%CV <sup>1</sup>	%背景 <sup>2</sup>
TG(56:6) [SIM]	924.8 → 924.8	TG(56:6) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.5	<1
TG(56:7) [NL-20:4]	922.8 → 601.5	TG(56:7) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.6	<1
TG(56:7) [NL-20:5]	922.8 → 603.5	TG(56:7) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.9	<1
TG(56:7) [NL-22:5]	922.8 → 575.5	TG(56:7) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.6	<1
TG(56:7) [NL-22:6]	922.8 → 577.5	TG(56:7) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	3.1	<1
TG(56:7) [SIM]	922.8 → 922.8	TG(56:7) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	2.2	<1
TG(56:8) [NL-20:4]	920.8 → 599.5	TG(56:8) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	3.2	<1
TG(56:8) [NL-20:5]	920.8 → 601.5	TG(56:8) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	3.7	<1
TG(56:8) [NL-22:6]	920.8 → 575.5	TG(56:8) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	2.4	<1
TG(56:8) [SIM]	920.8 → 920.8	TG(56:8) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	2.3	<1
TG(56:9) [NL-22:6]	918.8 → 573.5	TG(56:9) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	3.1	<1
TG(56:9) [SIM]	918.8 → 918.8	TG(56:9) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	2.6	57.34
TG(58:10) [NL-22:6]	944.9 → 599.5	TG(58:10) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	3.7	<1
TG(58:10) [SIM]	944.9 → 944.9	TG(58:10) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	2.2	<1
TG(58:8) [NL-22:6]	948.8 → 603.5	TG(58:8) [NL-22:6] [reference]	TG(15:0_18:1-d7_15:0)	1	3.8	<1
TG(58:9) [NL-22:6]	946.9 → 601.5	TG(58:9) [SIM] [reference]	TG(15:0_18:1-d7_15:0)	1	4.0	<1
TG(O-50:1) [NL-15:0]	836.8 → 577.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	6.0	<1
TG(O-50:1) [NL-16:0]	836.8 → 563.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.1	<1
TG(O-50:1) [NL-17:1]	836.8 → 551.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	7.8	<1
TG(O-50:1) [NL-18:1]	836.8 → 537.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.9	<1
TG(O-50:1) [SIM]	836.8 → 836.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.6	<1
TG(O-50:2) [NL-16:1]	834.8 → 563.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	5.5	12.07
TG(O-50:2) [NL-18:1]	834.8 → 535.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	4.6	<1
TG(O-50:2) [NL-18:2]	834.8 → 537.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	7.0	<1
TG(O-50:2) [SIM]	834.8 → 834.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.1	35.14
TG(O-52:0) [NL-16:0]	866.8 → 593.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	8.3	2.04
TG(O-52:1) [NL-16:0]	864.8 → 591.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.3	<1
TG(O-52:1) [NL-18:1]	864.8 → 565.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	2.7	<1
TG(O-52:1) [SIM]	864.8 → 864.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	2.5	<1
TG(O-52:2) [NL-16:0]	862.8 → 589.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	4.5	<1
TG(O-52:2) [NL-17:1]	862.8 → 577.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	5.6	<1
TG(O-52:2) [NL-18:1]	862.8 → 563.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	4.7	<1
TG(O-52:2) [SIM]	862.8 → 862.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.3	<1
TG(O-54:2) [NL-17:1]	890.8 → 605.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.7	<1
TG(O-54:2) [NL-18:1]	890.8 → 591.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.0	<1
TG(O-54:2) [SIM]	890.8 → 890.8	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	3.3	<1
TG(O-54:3) [NL-17:1]	888.8 → 603.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	7.4	<1
TG(O-54:3) [NL-18:1]	888.8 → 589.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	4.0	<1
TG(O-54:4) [NL-17:1]	886.8 → 601.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	11.7	<1
TG(O-54:4) [NL-18:2]	886.8 → 589.5	TG(48:1) [NL-18:1] d7 (IS) [reference]	TG(15:0_18:1-d7_15:0)	1	5.6	<1
Ubiquinone	880.7 → 197.0	Ubiquinone [reference]	TG(15:0_18:1-d7_15:0)	1	2.3	<1

<sup>1</sup> 通过依次进样的 50 个 TQCs 计算每种脂质浓度的 %CV

<sup>2</sup> % 背景是指在相同 LC/MS 批次中，空白浓度中值占 NIST1950 平均浓度的百分比

查找当地的安捷伦客户中心：

[www.agilent.com/chem/contactus-cn](http://www.agilent.com/chem/contactus-cn)

免费专线：

800-820-3278, 400-820-3278 (手机用户)

联系我们：

[LSCA-China\\_800@agilent.com](mailto:LSCA-China_800@agilent.com)

在线询价：

[www.agilent.com/chem/erfq-cn](http://www.agilent.com/chem/erfq-cn)

[www.agilent.com](http://www.agilent.com)

仅供科研使用。不用于临床诊断用途。

RA44413.1612962963

本文中的信息、说明和指标如有变更，恕不另行通知。

© 安捷伦科技（中国）有限公司，2021  
2021年9月23日，中国出版  
5994-3747ZHCN

