

KEY FINDING

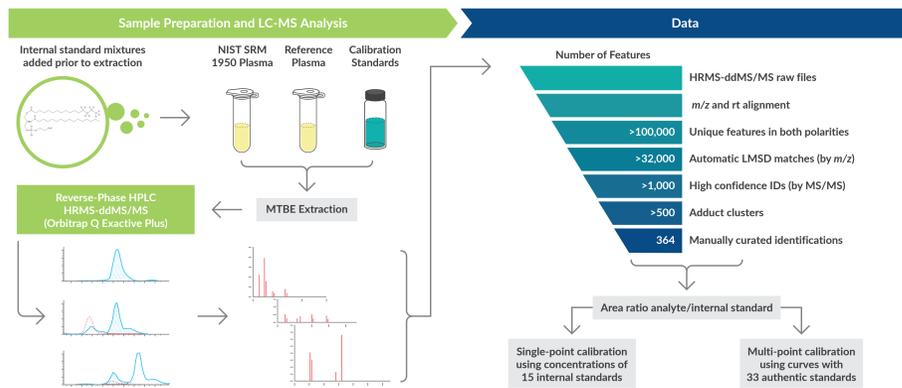
Untargeted lipidomics analysis provides useful quantitation of hundreds of molecular species in human plasma using single-point or multipoint calibration, the latter being more accurate across all lipid classes.

INTRODUCTION

- Untargeted lipidomics uses MS to analyze hundreds of molecular species in biological samples.
- Single-point calibration based on the known amounts of internal standards is used in many studies, especially those using shotgun (i.e., without chromatography) MS analysis.
- Questions remain about the accuracy of single-point calibration compared to the traditional interpolation in multipoint calibration curves, as well as its suitability in LC-MS studies, where internal standards may not coelute with most of the endogenous analytes.
- The objective of this study is to assess the accuracy of quantitation of lipids in human plasma by LC-MS using either single-point or multipoint calibration curves with authentic or surrogate lipid standards, using a new preparation of human reference plasma and methods validated by comparison with published values from the NIST SRM 1950 Metabolites in Human Plasma material.

(The authors declare no competing financial interest.)

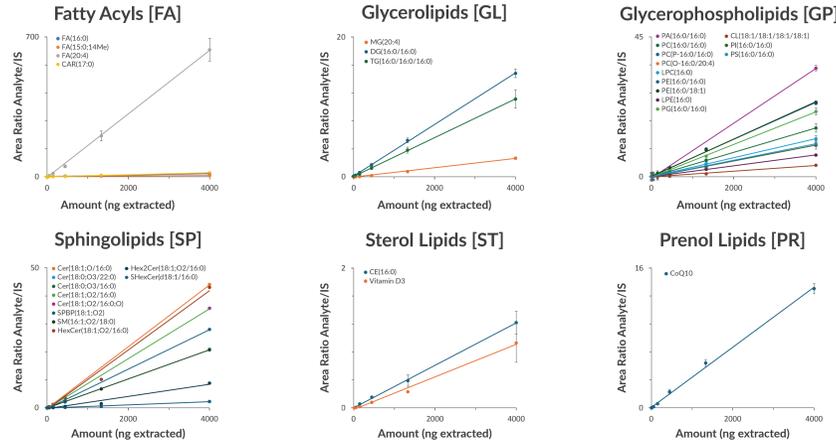
WORKFLOW



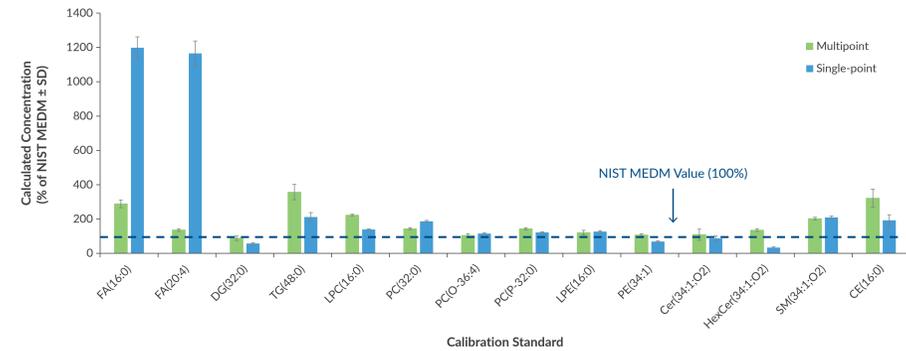
Calibration Standard	Internal Standard Used	LLOQ (ng)	ULOQ (ng)	Internal Standard	Amount Extracted (ng)
FA(16:0)	FA(16:0-d)	49.4	4000	FA(16:0-d)	400
FA(15:0,14Me)	FA(16:0-d)	49.4	4000	CAR(17:0-d)	50
FA(20:4(5Z,8Z,11Z,14Z))	FA(16:0-d)	1.82	4000	DG(16:0-d ₁ /16:0/0:0)	50
CAR(17:0)	CAR(17:0)-d ₅	0.2	4000	TG(16:0-d ₁ /16:0/16:0)	100
MG(20:4)	TG(16:0-d ₁ /16:0/16:0)	5.4	4000	PC(16:0-d ₁ /16:0)	100
DG(16:0/16:0)	DG(16:0-d ₁ /16:0)	0.6	4000	LPC(16:0-d ₁)	100
TG(16:0/16:0/16:0)	TG(16:0-d ₁ /16:0/16:0)	0.2	4000	PE(16:0-d ₁ /16:0)	50
PA(16:0/16:0)	PE(16:0-d ₁ /16:0)	5.4	4000	LPE(16:0-d ₁ /0:0)	100
PC(16:0/16:0)	PC(16:0-d ₁ /16:0)	0.6	4000	PG(16:0-d ₁ /16:0)	100
PC(P-16:0/16:0)	PC(16:0-d ₁ /16:0)	1.8	4000	PI(16:0-d ₁ /16:0)	100
PC(O-16:0/20:4)	PC(16:0-d ₁ /16:0)	5.4	4000	PS(16:0-d ₁ /16:0)	100
LPC(16:0)	LPC(16:0-d ₁)	0.6	4000	Cer(d18:1-d ₁ /16:0)	50
PE(16:0/16:0)	PE(16:0-d ₁ /16:0)	0.6	4000	SM(d18:1/16:0-d ₁)	100
PE(16:0/18:1(9Z))	PE(16:0-d ₁ /16:0)	0.6	4000	Cholesterol-d ₅	1000
LPE(16:0)	LPE(16:0-d ₁)	0.2	4000	CE(16:0-d ₁)	1000
PG(16:0/16:0)	PG(16:0-d ₁ /16:0)	0.2	4000		
PLG(18:0)	PE(16:0-d ₁ /16:0)	0.6	4000		
CL(18:1(9Z)/18:1(9Z)/18:1(9Z)/18:1(9Z))	PE(16:0-d ₁ /16:0)	5.4	4000		
PI(16:0/16:0)	PI(16:0-d ₁ /16:0)	0.6	4000		
PS(16:0/16:0)	PS(16:0-d ₁ /16:0)	5.4	4000		
Cer(18:1/O/16:0)	Cer(d18:1-d ₁ /16:0)	1.8	4000		
Cer(18:0/O3/22:0)	Cer(d18:1-d ₁ /16:0)	1.8	4000		
Cer(18:0/O3/16:0)	Cer(d18:1-d ₁ /16:0)	1.8	4000		
Cer(18:1/O2/16:0)	Cer(d18:1-d ₁ /16:0)	0.6	4000		
Cer(18:1/O2/16:0/O)	Cer(d18:1-d ₁ /16:0)	0.6	4000		
SPBP(18:1/O2)	Cer(d18:1-d ₁ /16:0)	50	4000		
SM(16:1;O2/18:0)	SM(d18:1/16:0-d ₁)	1.8	4000		
HexCer(18:1;O2/16:0)	Cer(d18:1-d ₁ /16:0)	0.6	4000		
Hex2Cer(18:1;O2/16:0)	Cer(d18:1-d ₁ /16:0)	1.8	4000		
SHexCer(18:1/16:0)	Cer(d18:1-d ₁ /16:0)	0.6	4000		
CE(16:0)	CE(16:0-d ₁)	49.4	4000		
Vitamin D3	TG(16:0-d ₁ /16:0/16:0)	49.4	4000		
CoQ10	TG(16:0-d ₁ /16:0/16:0)	0.6	4000		

RESULTS

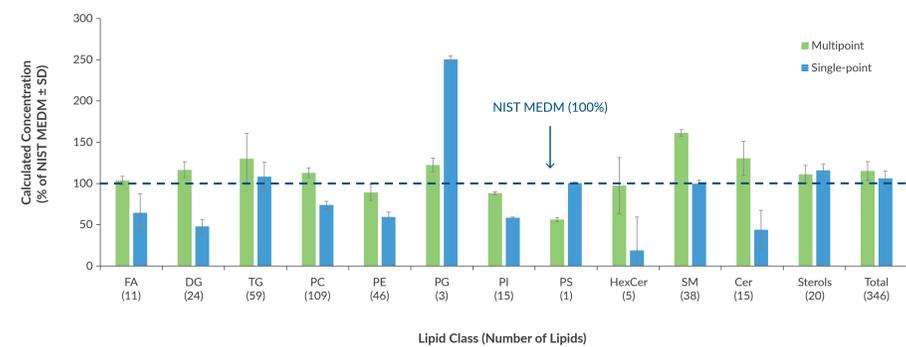
Calibration curves used for multipoint calculations



Validation of the calculated concentrations using single-point or multipoint calibration, as compared with the published median of the means (MEDM) for NIST SRM 1950 plasma



Comparison of total concentrations of each lipid class, as the sum of all molecular species identified, using single-point or multipoint calibration



Calculated concentrations (ng/μl) of the sums of glycerophosphocholines using single-point calibration or multipoint calibrations using the curves for three different PC standards

PC Subclass (number of features)	Published MEDM (ng/μl)	Curve Used for Multipoint Calibration			Single-Point Calibration
		PC(16:0/16:0)	PC(O-16:0/20:4)	PC(P-16:0/16:0)	
Diacylglycerophosphocholines (18)	502.9	502.9	502.9	502.9	686.0
1-Alkyl,2-aclyglycerophosphocholines (6)	31.6	502.9	502.9	502.9	33.8
1-(1Z-Alkenyl),2-aclyglycerophosphocholines (6)	20.7	502.9	502.9	502.9	18.0
Relative Error (%)	-	17.3	157.1	107.1	32.9

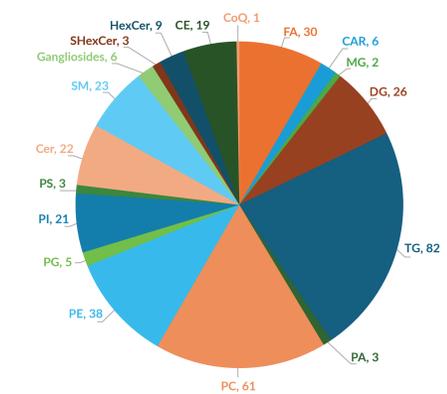


Download this poster and related information from our website

Effect of volume extracted in the number of curated molecular species quantitated

Lipid Class	Volume of Plasma Extracted (μL)				
	60	20	6.7	2.2	0.74
FA	30	30	27	24	6
CAR	6	6	2	0	0
MG	2	2	2	2	0
DG	28	26	26	27	14
TG	82	82	82	77	36
PA	3	3	3	1	0
PC	63	61	56	49	27
PE	38	38	36	26	11
PG	6	5	5	3	1
PI	21	21	20	18	9
PS	3	3	2	0	0
Cer	22	22	21	18	4
\$1P	1	0	0	0	0
SM	23	23	23	20	8
Gangliosides	6	6	6	6	0
SHexCer	3	3	3	2	0
HexCer	10	9	8	8	3
Sterols	21	19	15	15	6
CoQ10	1	1	1	0	0
Total	369	360	338	296	125

Number of curated molecular species identified by class



Total lipid concentration (ng/μl), by class

Lipid Class (Abbreviation)	Multipoint Calibration	
	Concentration	SD
Fatty Acids and Conjugates (FA)	380.4	43.2
Monoradylglycerols (MG)	0.5	0.0
Diradylglycerols (DG)	110.5	10.6
Triradylglycerols (TG)	647.0	107.0
Glycerophosphates (PA)	2.2	0.2
Glycerophosphocholines (PC)	804.9	20.2
Glycerophosphoethanolamines (PE)	24.2	2.1
Glycerophosphoglycerols (PG)	15.7	1.1
Glycerophosphoinositols (PI)	46.5	0.2
Glycerophosphoserines (PS)	1.0	0.0
Ceramides (Cer)	14.6	1.2
Phosphosphingolipids (SM)	420.5	20.5
Gangliosides	0.2	0.0
Sulfatides (SHexCer)	0.1	0.0
Neutral glycosphingolipids (HexCer)	6.4	1.0
Sterols (CE)	5399.0	175.6
Quinones (CoQ)	0.8	0.1
Total	7874.4	381.0

CONCLUSIONS

- Untargeted lipidomics analysis by LC-MS provides excellent linearity of the signals for 33 lipid standards across several orders of magnitude.
- Calculated plasma concentrations using either single-point or multipoint calibration are in reasonable agreement with published data, validating the workflow.
- Although there are differences dependent on the lipid classes monitored, multipoint calibration results in improved accuracy across all the classes.
- Better results are obtained when using the isotopically labeled internal standard corresponding to the calibration standard, and a calibration curve from a structurally similar standard.
- The method presented here allows for the confident identification and quantitation of hundreds of lipid molecular species from sub-microliter volumes of human plasma.