

Application Note

Characterization of Human Plasma as a Reference Material for Lipid Analysis Using LC-MS

Matthew J. Kwiatkowski, Samantha K. Goodwin, Samuel L. DeLoy, Dilrukshika S.W. Palagama, Miguel A. Gijón, and Paul D. Kennedy
Cayman Chemical, Ann Arbor, MI

Key Features

- A preparation of human plasma has been characterized as a reference material for lipid identification and quantitation using LC-MS.
- The validity of the methods used and the values obtained is supported by the reasonable agreement with published values in NIST SRM 1950 plasma.
- MaxSpec® Reference Plasma (Human) can be used for quality control, method development, and performance validation in lipid studies.
- A list of lipid concentrations is made public and periodically updated.

To cite this application note:

Kwiatkowski, M.J., Goodwin, S.K., DeLoy, S.L., *et al.* Characterization of human plasma as a reference material for lipid analysis using LC-MS. *Application Note, Cayman Chemical Company* (2025).

Introduction

Lipids are key molecules for maintaining cellular membranes, energy storage, cellular signaling, and many other biological processes. The identification and understanding of their diverse functions are useful for insights into cell biology, metabolism, and disease mechanisms. Mass spectrometry has become the standard analytical tool for the identification and quantitation of hundreds of lipids, helping to reveal their roles in biological systems.

The use of well-characterized reference materials increases confidence in lipid quantitative data by validating the analytical methods employed. The objective of this study is to characterize a preparation of human plasma so it can be made available as reference material for the analysis of a wide variety of lipids using targeted methods for the analysis of oxylipins and bile acids, as well as untargeted lipidomics methods. All methods used in this characterization are validated by comparison of the obtained values using the well-established NIST SRM 1950 Metabolites in Human Plasma material to the published values.^{1,2}

Experimental Procedures



Figure 1. Workflow of lipid extraction and analysis of plasma using LC-MS.

Replicate 50 µl aliquots (three for each analytical method) of each plasma (MaxSpec® or NIST SRM 1950) were extracted after mixtures of deuterated internal standards were added (See Table 1 for internal standard list). Acetonitrile extraction was used to precipitate proteins and isolate bile acids.³ Oxylipins were isolated using reversed-phase polymeric sorbent solid-phase extraction (SPE).⁴ For untargeted analysis, methyl-tert-butyl ether-based extraction was used.⁵ Separation and quantitation of bile acids and oxylipins were achieved by reversed-phase LC-MS/MS using an Exion UPLC system coupled with a triple-quadrupole 6500+ mass spectrometer (Sciex). Separation and quantitation of all lipids using an untargeted approach were achieved using reversed-phase HPLC coupled with a high-resolution Q Exactive™ Orbitrap™ mass spectrometer (Thermo Fisher Scientific).

Table 1. Deuterated internal standards used for the three types of analysis. ¹Components of Deuterated Bile Acids MaxSpec® Discovery Mixture; Item No. 33506. ²Components of Deuterated Lipidomics MaxSpec® Mixture; Item No. 40974.

| Oxylipins | Item No. | Bile Acids ¹ | Item No. | Untargeted ² | Item No. |
|---|----------|---|----------|--|----------|
| PGE ₂ -d ₄ | 10007273 | Cholic Acid-d ₄ | 31348 | FA(16:0-d ₅) | 30557 |
| TXB ₂ -d ₄ | 319030 | Taurocholic Acid-d ₄ | 31375 | CAR(17:0)-d ₃ | 9003263 |
| 5-HETE-d ₈ | 334230 | Glycocholic Acid-d ₄ | 31352 | DG(16:0-d ₉ /16:0) | 27591 |
| LTB ₄ -d ₄ | 320110 | Deoxycholic Acid-d ₄ | 31350 | TG(16:0-d ₉ /16:0/16:0) | 27592 |
| LTD ₄ -d ₅ | 10006199 | Taurodeoxycholic Acid-d ₄ | 31563 | PC(16:0-d ₉ /16:0) | 28154 |
| LXA ₄ -d ₅ | 10007737 | Glycodeoxycholic Acid-d ₄ | 31553 | LPC(16:0-d ₉) | 28153 |
| 14(15)-EET-d ₁₁ | 10006410 | Chenodeoxycholic Acid-d ₄ | 31366 | PE(16:0-d ₉ /16:0) | 28155 |
| 14(15)-DiHET-d ₁₁ | 10008040 | Taurochenodeoxycholic Acid-d ₄ | 31362 | LPE(16:0-d ₉) | 27588 |
| 9-HODE-d ₄ | 338410 | Glycochenodeoxycholic Acid-d ₄ | 31364 | PG(16:0-d ₉ /16:0) | 28953 |
| RvE1-d ₄ | 10009854 | Lithocholic Acid-d ₄ | 31354 | PI(16:0-d ₉ /16:0) | 28156 |
| RvD1-d ₅ | 11182 | Taurolithocholic Acid-d ₄ | 31571 | PS(16:0-d ₉ /16:0) | 28152 |
| 12,13-DiHOME-d ₄ | 10009994 | Glycolithocholic Acid-d ₄ | 31554 | Cer(18:1(4E)-d ₇ ;1OH,3OH/16:0) | 22787 |
| 13,14-Dihydro-15kPGD ₂ -d ₉ | 19334 | Ursodeoxycholic Acid-d ₄ | 31368 | SM(18:1(4E)-;1OH,3OH/16:0-d ₉) | 27551 |
| 15-HETE-d ₈ | 334720 | Tauroursodeoxycholic Acid-d ₄ | 31564 | Cholesterol-d ₇ | 25265 |
| PGA ₂ -d ₄ | 310210 | Glycoursodeoxycholic Acid-d ₄ | 31555 | CE(16:0-d ₉) | 28123 |

To quantitate bile acids and oxylipins, the integrated area of the LC-MS/MS signal corresponding to each analyte was normalized to the integrated area of its corresponding internal standard, when possible, or a surrogate internal standard. Using MultiQuant software (Sciex), these area ratios were interpolated into an authentic or surrogate calibration curve to provide a calculated concentration value. Untargeted lipidomics data was analyzed using Lipostar software for feature detection, peak alignment, normalization, and lipid identification. The integrated areas corresponding to all detected analytes were also normalized to the integrated areas corresponding to internal standards within their class, when possible, and single-point calibration using the known amount of internal standard added to samples was used to calculate the concentrations of all analytes. In this initial study, only those lipid molecular species for which published data were available for validation using the NIST plasma were quantified.

Results

Analysis of the NIST SRM 1950 plasma and a comparison of the data obtained with published NIST SRM values is a necessary initial step to validate the methods used to characterize the MaxSpec® plasma.^{1,2} As shown in Figure 2 and Table 2, the lipid analytical data compare well with published reference values for the NIST plasma.

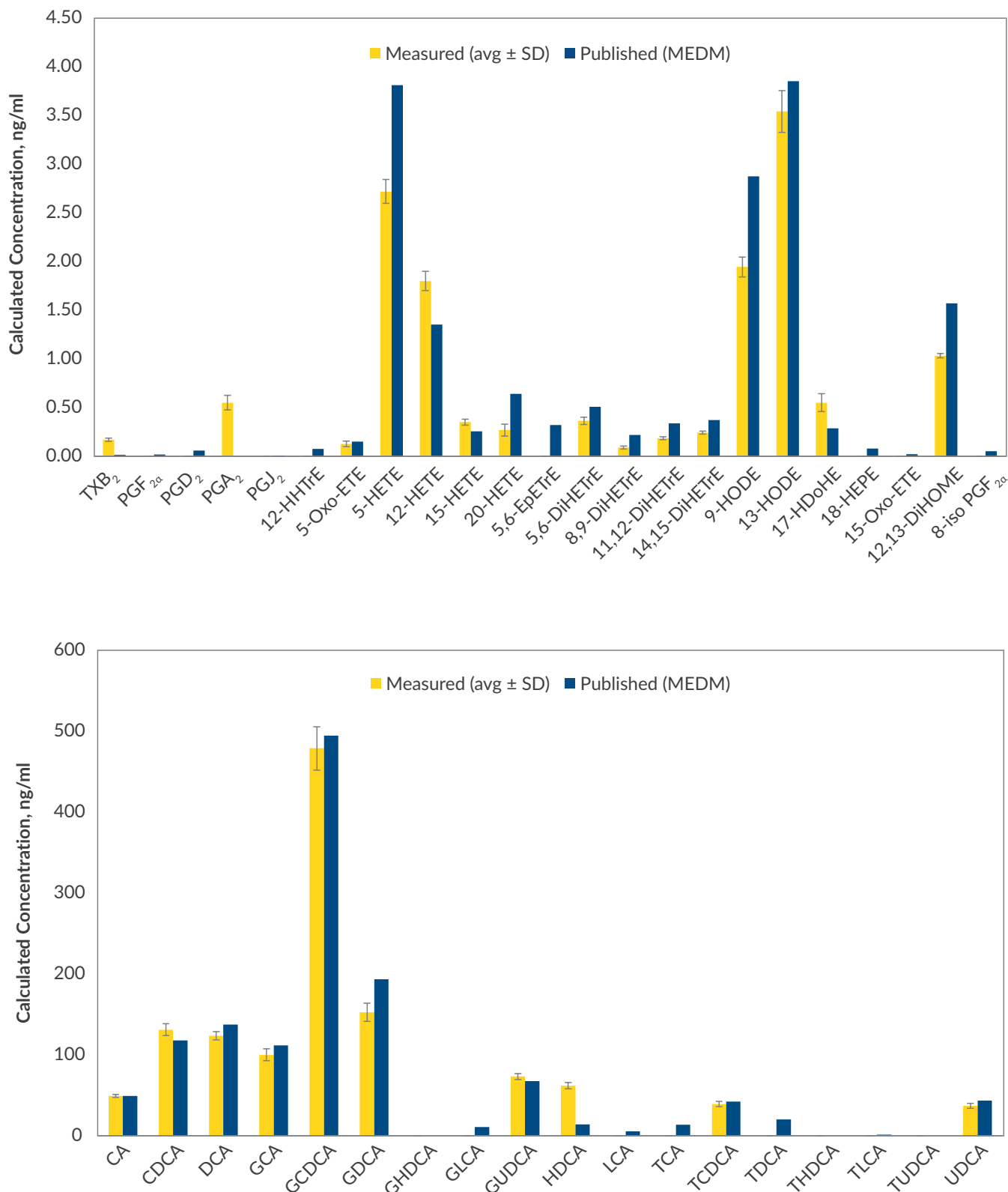


Figure 2. Validation of the methods used for targeted analysis of oxylipins (top) and bile acids (bottom) in NIST reference plasma SRM 1950. The measured data is presented as avg ± SD (n=15; 3 replicate extractions, 5 injections each) and the published data is the consensus MEDM (median of the means).

Table 2. Validation of the method used for untargeted lipidomics analysis in NIST reference plasma SRM 1950. Data shown are the percent of each molecular species in its class, with only the five most abundant lipid molecular species represented for clarity. Values shown are the averages of 12 replicates (3 replicate extractions, 4 injections each).

| Molecular Species (Sum Composition) | Published Values (MEDM) | Measured Values (This Study) |
|-------------------------------------|-------------------------|------------------------------|
| CAR(18:1) | N/A | 37.9 |
| CAR(18:2) | N/A | 18.5 |
| CAR(16:0) | N/A | 17.9 |
| CAR(14:1) | N/A | 9.4 |
| CAR(18:0) | N/A | 6.9 |
| DG(36:3) | 15.7 | 18.6 |
| DG(34:0) | 12.2 | 6.6 |
| DG(36:2) | 11.6 | 13.4 |
| DG(34:1) | 11.4 | 18.8 |
| DG(34:2) | 8.2 | 14.9 |
| TG(52:3) | 16.9 | 15.1 |
| TG(52:4) | 8.1 | 10.6 |
| TG(50:2) | 7.9 | 7.5 |
| TG(52:2) | 7.4 | 14.0 |
| TG(50:1) | 6.4 | 5.3 |
| LPC(16:0) | 46.4 | 43.9 |
| LPC(18:0) | 17.2 | 13.5 |
| LPC(18:2) | 14.0 | 12.6 |
| LPC(18:1) | 11.4 | 10.4 |
| LPC(20:4) | 3.8 | 2.9 |
| PC(34:2) | 20.3 | 34.9 |
| PC(36:4) | 12.7 | 1.6 |
| PC(36:2) | 11.8 | 22.7 |
| PC(34:1) | 10.2 | 2.2 |
| PC(36:3) | 8.5 | 13.9 |

| Molecular Species (Sum Composition) | Published Values (MEDM) | Measured Values (This Study) |
|-------------------------------------|-------------------------|------------------------------|
| LPE(18:2) | 23.8 | 29.4 |
| LPE(18:0) | 20.0 | 21.8 |
| LPE(18:1) | 17.5 | 15.8 |
| LPE(20:4) | 13.8 | 9.8 |
| LPE(16:0) | 11.4 | 13.1 |
| PE(38:4) | 15.1 | 7.5 |
| PE(36:2) | 12.5 | 6.0 |
| PE(O-38:7) | 6.5 | 2.7 |
| PE(38:6) | 6.0 | 7.1 |
| PE(36:4) | 5.8 | 2.0 |
| PI(38:4) | 39.5 | 34.1 |
| PI(36:2) | 16.0 | 23.2 |
| PI(38:3) | 7.1 | 8.7 |
| PI(36:4) | 6.2 | 6.4 |
| PI(34:2) | 5.8 | 6.2 |
| Cer(42:1;O2) | 34.9 | 39.3 |
| Cer(42:2;O2) | 15.1 | 19.0 |
| Cer(41:1;O2) | 12.3 | 11.4 |
| Cer(40:1;O2) | 12.0 | 12.6 |
| Cer(34:1;O2) | 5.1 | 3.4 |
| SM(34:1;O2) | 30.0 | 31.5 |
| SM(42:2;O2) | 13.2 | 14.5 |
| SM(36:1;O2) | 6.0 | 5.7 |
| SM(40:1;O2) | 6.0 | 8.5 |
| SM(42:1;O2) | 6.0 | 5.7 |
| CE(18:2) | 55.4 | 49.6 |
| CE(18:1) | 14.7 | 9.5 |
| CE(20:4) | 11.4 | 25.8 |
| CE(16:0) | 6.8 | 1.2 |
| CE(16:1) | 3.3 | 1.0 |

The three validated methods were utilized to characterize the MaxSpec® reference plasma. Tables 3 and 4 show the calculated concentrations of bile acids and oxylipins detected in this reference plasma, whereas Figure 3 summarizes the distribution of hundreds of molecular species of lipids in this material.

A list of quantified lipid molecular species in this study is available [online](#).

Table 3. Calculated concentrations of bile acids in MaxSpec® reference plasma. (B)LLOQ, (below) lower limit of quantitation.

| Bile Acids | Shorthand | Calibration Standard (Item No.) | LLOQ (ng/ml) | Concentration (ng/ml) Avg ± SD (n=15) |
|----------------------------|-----------|---------------------------------|--------------|--|
| Cholic Acid | CA | 31347 | 3.8 | 108.6 ± 7.58 |
| Chenodeoxycholic Acid | CDCA | 31365 | 20.8 | 235.2 ± 15.15 |
| Deoxycholic Acid | DCA | 31349 | 14.1 | 189.7 ± 7.8 |
| Glycocholic Acid | GCA | 31351 | 8.4 | 28.3 ± 1.38 |
| Glycochenodeoxycholic Acid | GCDCA | 31363 | 7.7 | 173.3 ± 7.55 |
| Glycodeoxycholic Acid | GDCA | 31599 | 7.4 | 58.1 ± 4.52 |
| Glycohyodeoxycholic Acid | GHDCA | 31600 | 14.1 | BLLOQ |
| Glycolithocholic Acid | GLCA | 31601 | 14.9 | BLLOQ |
| Glycoursodeoxycholic Acid | GUDCA | 31602 | 8.1 | 14.8 ± 1.46 |
| Hyodeoxycholic Acid | HDCA | 31606 | 40.5 | 40.7 ± 5.31 |
| Lithocholic Acid | LCA | 31353 | 15.3 | BLLOQ |
| Taurocholic Acid | TCA | 31374 | 33.8 | BLLOQ |
| Taurochenodeoxycholic Acid | TCDCA | 31361 | 31.2 | BLLOQ |
| Taurodeoxycholic Acid | TDCA | 31603 | 31.2 | BLLOQ |
| Taurohyodeoxycholic Acid | THDCA | 31614 | 1.9 | BLLOQ |
| Taurolithocholic Acid | TLCA | 31604 | 60.5 | BLLOQ |
| Tauroursodeoxycholic Acid | TUDCA | 31605 | 67 | BLLOQ |
| Ursodeoxycholic Acid | UDCA | 31367 | 14.1 | 17.6 ± 1.8 |

Table 4. Calculated concentrations of oxylipins in MaxSpec® reference plasma. (B)LLOQ, (below) lower limit of quantitation; *Surrogate standard.

| Oxylipins | Shorthand | Calibration Standard (Item No.) | LLOQ (ng/ml) | Concentration (ng/ml) Avg ± SD (n=15) |
|--|-------------------------|---------------------------------|--------------|--|
| Thromboxane B ₂ | TXB ₂ | 19030 | 0.04 | 3.32 ± 0.17 |
| Prostaglandin F _{2α} | PGF _{2α} | 10007211* | 0.12 | 1 ± 0.1 |
| Prostaglandin D ₂ | PGD ₂ | 10007211* | 0.12 | 0.41 ± 0.04 |
| Prostaglandin A ₂ | PGA ₂ | 18500* | 0.12 | 24.49 ± 0.97 |
| Prostaglandin J ₂ | PGJ ₂ | 18500 | 0.12 | 3.17 ± 0.23 |
| 12-Hydroxy-heptadecatrienoic Acid | 12-HHTrE | 34590 | 0.12 | 1.03 ± 0.08 |
| 5-Oxo-eicosatetraenoic Acid | 5-Oxo-ETE | 34250 | 0.04 | 132.49 ± 6.19 |
| 5-Hydroxy-eicosatetraenoic Acid | 5-HETE | 34210 | 0.01 | 2738 ± 101.11 |
| 12-Hydroxy-eicosatetraenoic Acid | 12-HETE | 34550 | 0.04 | 1477 ± 105.25 |
| 15-Hydroxy-eicosatetraenoic Acid | 15-HETE | 34550* | 0.04 | 1126 ± 59.57 |
| 20-Hydroxy-eicosatetraenoic Acid | 20-HETE | 10007269 | 0.04 | 0.71 ± 0.13 |
| 5,6-Epoxy-eicosatrienoic Acid | 5,6-EpETrE | 50211 | 0.37 | BLLOQ |
| 8,9-Epoxy-eicosatrienoic Acid | 8,9-EpETrE | 50211* | 0.37 | BLLOQ |
| 11,12-Epoxy-eicosatrienoic Acid | 11,12-EpETrE | 50511 | 0.01 | BLLOQ |
| 5,6-Dihydroxy-eicosatrienoic Acid | 5,6-DiHETrE | 51211 | 0.01 | 2.44 ± 0.09 |
| 8,9-Dihydroxy-eicosatrienoic Acid | 8,9-DiHETrE | 51211* | 0.01 | 2.88 ± 0.1 |
| 11,12-Dihydroxy-eicosatrienoic Acid | 11,12-DiHETrE | 51511 | 0.01 | 1.01 ± 0.05 |
| 14,15-Dihydroxy-eicosatrienoic Acid | 14,15-DiHETrE | 51511* | 0.01 | 2.12 ± 0.11 |
| 9-Hydroxy-octadecadienoic Acid | 9-HODE | 38400 | 0.01 | 2743 ± 139.27 |
| 13-Hydroxy-octadecadienoic Acid | 13-HODE | 38400* | 0.01 | 3213 ± 118.34 |
| 17-Hydroxy-docosahexaenoic Acid | 17-HDoHE | 33650 | 0.12 | 520 ± 31.56 |
| 18-Hydroxy-eicosapentaenoic Acid | 18-HEPE | 32840 | 0.04 | 22.46 ± 1.07 |
| 15-Oxo-eicosatetraenoic Acid | 15-Oxo-ETE | 34550* | 0.04 | 652.40 ± 38.71 |
| 12,13-Dihydroxy-octadecamonoenoic Acid | 12,13-DiHOME | 10009832 | 0.01 | 107.11 ± 3.24 |
| 9,10-Dihydroxy-octadecamonoenoic Acid | 9,10-DiHOME | 10009832* | 0.01 | 137.85 ± 4.36 |
| 8-Iso-prostaglandin F _{2α} | 8-iso-PGF _{2α} | 18500* | 0.12 | 2.5 ± 0.46 |

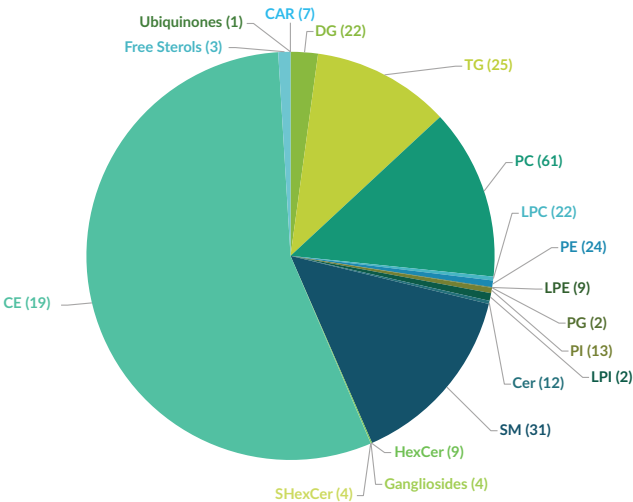


Figure 3. Distribution of lipid molecular species by their measured abundance in MaxSpec® reference plasma.

As additional characterization of both the MaxSpec® plasma and the untargeted lipidomics method, we explored the effect that decreasing volumes of plasma extracted would have in the number of molecular species identified. As expected, the total number of detectable molecular species is higher with higher volumes extracted, as shown in Table 5. However, the increase of identifications is relatively modest above 2 µl. Of note, the untargeted analysis utilized allows for the identification and quantitation of over a hundred lipid molecular species from sub-microliter volumes of human plasma.

Table 5. Effect of volume extracted on the number of curated molecular species detectable in the MaxSpec® reference plasma.

| Lipid Class | Volume of Plasma Extracted (µl) | | | | |
|--------------|---------------------------------|------------|------------|------------|------------|
| | 60 | 20 | 6.7 | 2.2 | 0.74 |
| 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 |
| S1P | 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 | 339 | 330 | 311 | 272 | 119 |

Additional characterization of this reference plasma preparation is ongoing to increase the coverage of quantified lipids and to evaluate the stability of analytes, which will progressively increase the usefulness of this material as a reference control.

References

1. Bowden, J.A., Heckert, A., Ulmer, C.Z., *et al.* Harmonizing lipidomics: NIST interlaboratory comparison exercise for lipidomics using SRM 1950-metabolites in frozen human plasma. *J. Lipid Res.* **58(12)**, 2275-2288 (2017).
2. Quehenberger, O., Armando, A.M., Brown, A.H., *et al.* Lipidomics reveals a remarkable diversity of lipids in human plasma. *J. Lipid Res.* **51(11)**, 3299-3305 (2010).
3. Krautbauer, S. and Liebisch, G. LC-MS/MS analysis of bile acids. *Methods Mol. Biol.* **1730**, 103-110 (2018).
4. Wang, Y., Armando, A.M., Quehenberger, O., *et al.* Comprehensive ultra-performance liquid chromatographic separation and mass spectrometric analysis of eicosanoid metabolites in human samples. *J. Chromatogr. A* **1359**, 60-69 (2014).
5. Lange, M., Angelidou, G., Ni, Z., *et al.* *AdipoAtlas*: A reference lipidome for human white adipose tissue. *Cell Rep. Med.* **2(10)**, 100407 (2021).