

(The authors declare no competing financial interest.)

- The targeted and untargeted quantitation methods used in this study produce robust lipid analytical data, as validated by the reasonable agreement with published reference values.
- The commercially-available reference plasma encompasses a broad spectrum of lipid molecular species and can serve as a valuable resource for validating lipid analysis methods.
- The reference plasma preparation contains surprisingly high amounts of several fatty acids, especially oxidized fatty acids. The reason for this is currently unknown but may be related to plasma collection or handling.
- 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 increase the usefulness of this material as a reference control.