

<sup>4</sup>Thermo Fisher Scientific, San Jose, CA, USA

- New nominal mass instruments with high acquisition rates, sensitivity, and advanced features like MSn are emerging.
- Proposal of an alternative to untargeted analysis on high-resolution instruments.
- Spectra will contain all product ions.
- PRM Conductor helps with product ion selection for data analysis.
- Large sex differences between male and female zebrafish livers indicate the importance of separating them in statistical data analysis.

- Targeted workflows using triple quadrupole (QQQ) instruments have recently gained more attention in the field of metabolomics [1].
- The targeted aspect, namely having to decide which metabolites to measure before data acquisition, seems to be a barrier for untargeted metabolomics
- New nominal mass instrument, Thermo Scientific™ Stellar™ Mass Spectrometer provides:
- Detection of all product ions, thus only precursors need to be known (can be identified or unknown compounds).
- High acquisition rates, and sensitivity.
- Features such as MSn, HCD, and CID, stepped collision energies.
- Software to help with method development.
- Zebrafish (*Danio rerio*) provides a premier model organism to study whole-body metabolism [2]. Using a targeted discovery metabolomics experiment, we measured liver extracts from adult male and female zebrafish.

# Sample preparation

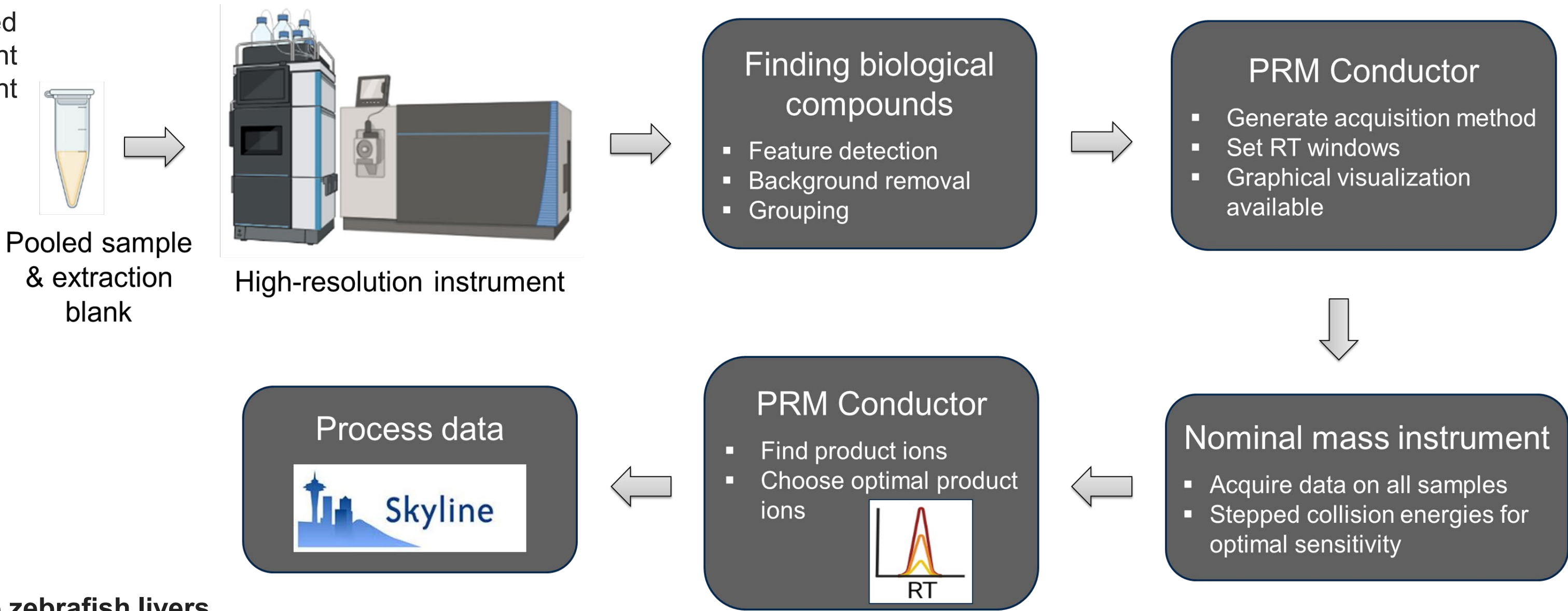
The flowchart illustrates the sample preparation process for lipid LC-MS analysis, starting from an anesthetized fish and ending with lipid LC-MS analysis. The steps are as follows:

- Anesthetize fish**: A small fish is shown being anesthetized.
- Collect individual tissues into pre-washed tubes**: Tissues are collected into a tube.
- Snap-freeze individual tissue in liquid N<sub>2</sub> and weigh**: The tissue is snap-frozen in liquid nitrogen and weighed.
- Grind with polypropylene pestle**: The tissue is ground using a pestle.
- Add 2.2 L mix according to recipe**: A liquid mixture is added to the ground tissue.
- 3x freeze-thaw, vortex 20" C x 3**: The mixture undergoes three freeze-thaw cycles and is vortexed at 20°C for 3 minutes.
- Centrifuge 14,000 g, 10 min, 4°C**: The mixture is centrifuged at 14,000 g for 10 minutes at 4°C.
- Supernatant**: The supernatant is collected from the centrifugation step.
- Polar LC-MS analysis**: The supernatant is analyzed using polar LC-MS.
- Re-extract pellet with 100% IPA**: The pellet is re-extracted with 100% isopropanol (IPA).
- 2x sonicate, vortex 20" C x 3**: The re-extracted pellet is sonicated twice and vortexed at 20°C for 3 minutes.
- Centrifuge 14,000 g, 10 min, 4°C**: The mixture is centrifuged again at 14,000 g for 10 minutes at 4°C.
- Mix 1:1 with pellet extract**: The supernatant from the second centrifugation is mixed 1:1 with the pellet extract.
- Lipid LC-MS analysis**: The final mixture is analyzed using lipid LC-MS.

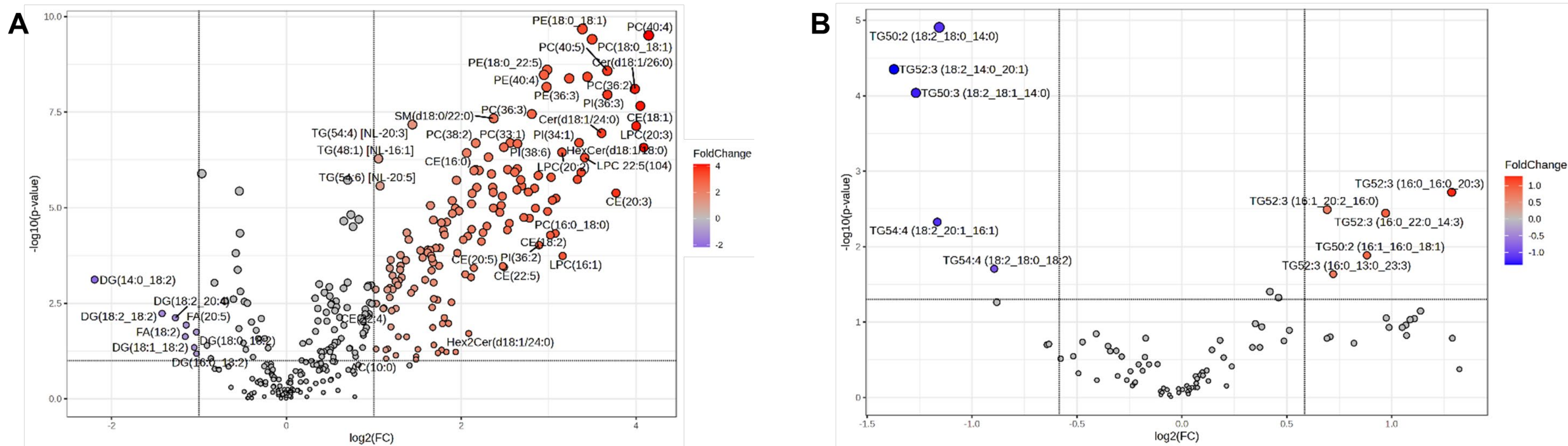
**Figure 1.** Sample preparation workflow for polar and lipid metabolite analysis from zebrafish.

- Peak list obtained from high-resolution instrument in both positive and negative polarity using Thermo Scientific™ Compound Discoverer™ 3.4 software.
- Removal of in-source fragments, filtering to keep only one polarity.
- List can contain identified and unidentified compounds.
- Thermo Scientific™ AcquireX in combination with Compound Discoverer 3.4 software and Thermo Scientific™ LipidSearch™ software used for the identification.
- Additional MS3 data can be acquired for triglycerides (TGs).

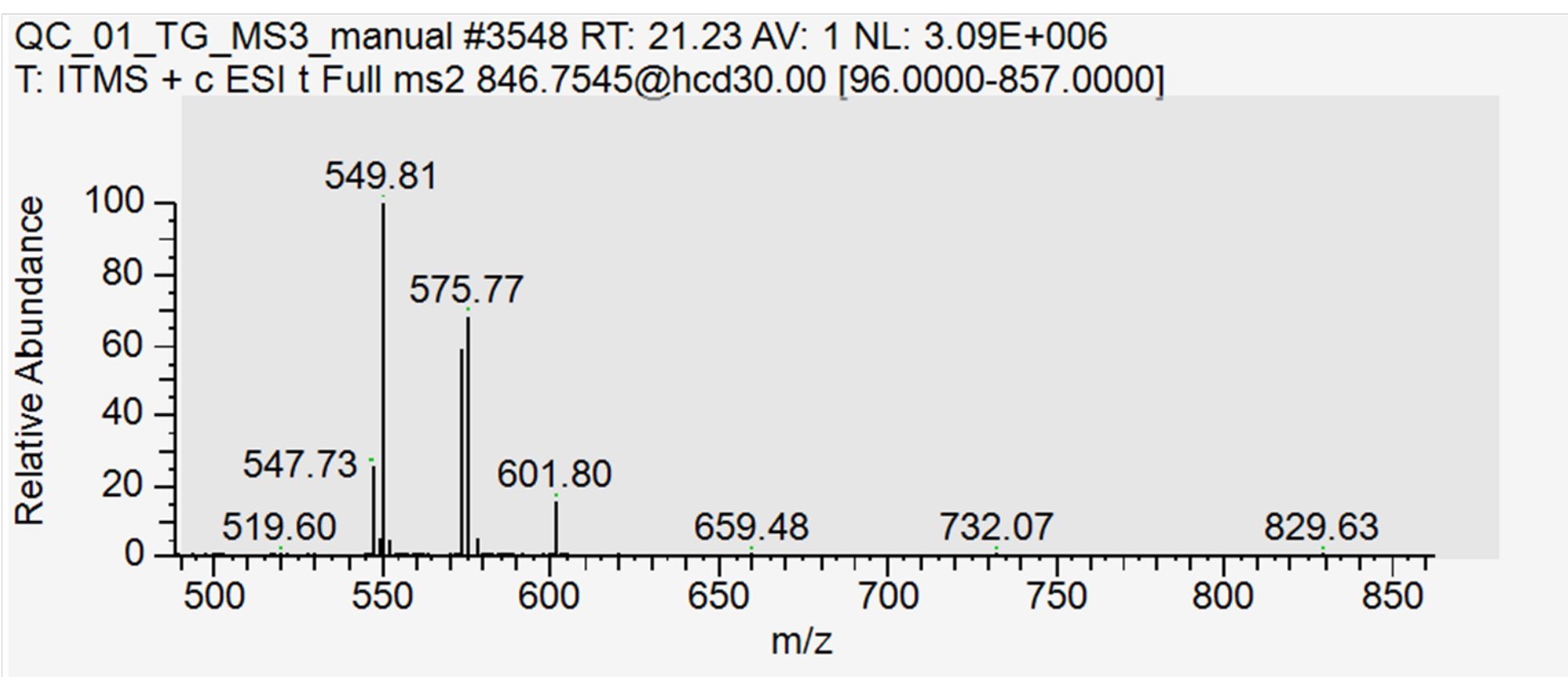
- Analysis of pooled sample and extraction blank on the high-resolution instrument
- Feature detection, grouping, etc. using Compound Discoverer 3.4 software
- PRM Conductor helps generate targeted acquisition methods with optimal RT windows
- Data acquisition for all samples on nominal mass instrument Stellar MS
- Stepped collision energies for optimal sensitivity, especially for unknowns
- Optimized collision energies can be used if available
- HCD and CID available (per precursor)
- PRM conductor helps to find optimal product ions and to generate a Skyline transition list
- Which product ions are the best (intensity and no interferences)



- Major differences in lipid metabolite levels between male and female livers.
- The majority of compounds were higher in females, but eight lipids were lower in females, seven of which contained FA 18:2.
- TGs containing FA 18:2 follow the same pattern (lower in females).
- TGs containing FA 18:2 and FA 14:0 show biggest difference.



**Figure 3.** Volcano plot of lipids detected in female vs. male zebrafish livers. A) lipids from MS2 analysis, B) MS3 analysis of selected TGs. Generated using MetaboAnalyst 6.0.

MS2 of TG(50:3) as  $[M+NH_4]^+$ 

**Figure 4.** MS2 and MS3 spectra of TG(50:3) as an example for triglyceride structural characterization.

- Streamlined approach from untargeted metabolomics to a targeted discovery methodology
- All product ions are detected, which can help distinguish isomers
- Large biological sex differences indicate the importance of analyzing data from each gender separately

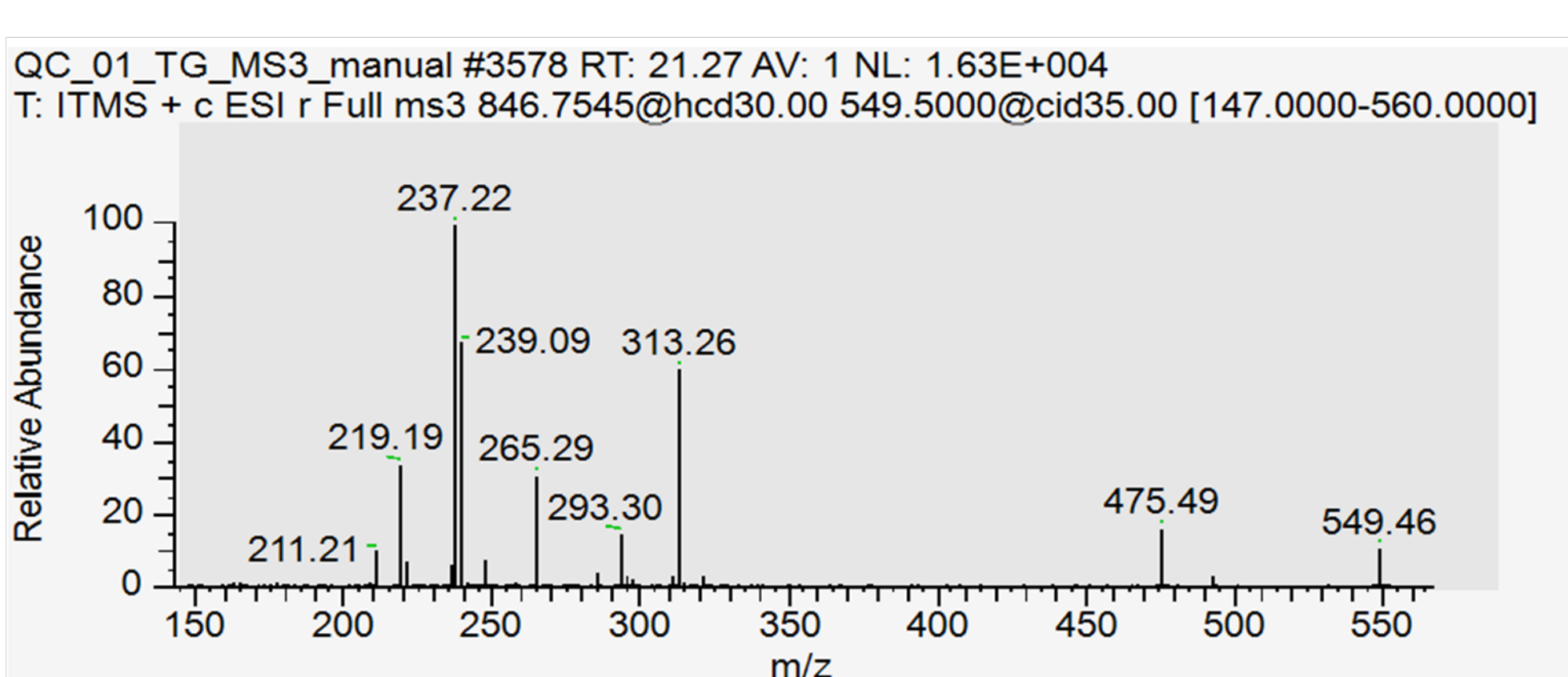
The authors thank Madelyn M. Jackstadt and Madison Barr for technical assistance with the zebrafish sample collection. This work was supported by funding from the National Institutes of Health grants U01CA235482 (G.J.P.) and R35ES028365 (G.J.P.).

Figure 2 was created using BioRender.com.

[1] Schwaiger-Haber et al., ACS Meas. Sci. 1, 1 (2021)

[2] Naser, Jackstadt et al., Cell Metab. 33, 7 (2021)

MS3 of 549



G.J.P. is a scientific advisory board member for Cambridge Isotope Laboratories and has a collaborative research agreement with Thermo Scientific. B.A., C.C.J., P.R., and S.S.B. work at Thermo Scientific. PO003608

All other authors do not have any conflicts of interest.

© 2025 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is not intended to encourage the use of these products in any manner that might infringe the intellectual property rights of others.