

Poster Reprint

ASMS 2018 MP-486

Informatics Solutions for Improved Identifications in 2D-LC/MS and IM-MS Lipidomics Workflows

Sarah M. Stow¹, Jeremy P. Koemel², Sonia Liggi³, Christine Hinz³, Julian L. Griffin³, Alex Apffel¹, Mark Sartain¹, Xiangdong Li¹, Norton Kitagawa¹, and John C. Fjeldsted¹

¹Agilent Technologies, Santa Clara, CA

²Department of Pathology, Immunology and Laboratory Medicine, University of Florida, Gainesville, FL

³Department of Biochemistry and Cambridge Systems Biology Centre, University of Cambridge, Cambridge, UK

Introduction

Lipidomics Workflows

Confident identifications are a challenge in the field of MS based lipidomics due to high occurrence of lipid isomers and isobars. Chromatographic separations prior to MS analysis can remove some signal overlap by separating lipid classes according to head group as well as separating lipids based on fatty acid saturation. Combining these two separations with a 2D-LC platform removes ambiguity for many ion signals. Performing data-independent MS/MS experiments after ion mobility (IM) separations allows fragment ions to be linked back to their precursor ion in the drift time dimension. These technologies offer key advances in confidently identifying more lipid species in omics workflows. In this study we apply a novel twist on the contemporary online high resolution 2D-LC experiment as well as evaluate the ability to make confident identifications with the All Ions IM-MS workflow with both 1D and 2D LC separations.

Two Dimensional Liquid Chromatography

1D and 2D LC experiments were performed on a commercial LC (1290 Series), Agilent Technologies, Santa Clara, CA. Multiple heart cutting and high resolution 2D LC experiments were performed with orthogonal LC methods (HILIC and Reverse Phase). The instrument was equipped with the 2 pos/4 port duo 2x6 loops.



Ion mobility mass spectrometry

The 6560 IM-Q-TOF (Agilent Technologies, Santa Clara, CA) was used for IM experiments. Single field CCS values were calculated for the LC-IM-MS experiments. Additionally, All Ions IM-MS fragmentation experiments were performed which align fragment ions with their precursors according to drift time.



LC Experiments

LC methods were adapted from a previously described method¹, but briefly, all first dimensional experiments were performed with an RX-Sil HILIC column (3.0 x 100 mm, 1.8 micron, 360 µL/min flow rate). Second dimension experiments were performed with a Zorbax SB-C18 column (2.1 x 100 mm, 1.8 micron, 0.2 mL/min flow rate). Ion mobility experiments were performed with the single field approach described previously². Lipid standards were purchased from Avanti Polar Lipids (Alabaster, AL) and were run as both a mixture of standards as well as spiked into a bovine heart extract (Avanti).

1D-LC			
Mi n	Sol A: ACN (0.1% FA)	Sol B: ACN:MeOH:H₂O (50:20:30 v/v) (20 mM NH₄HCO₂)	
0	70%	30%	
2	40%	60%	
4	30%	70%	
5	0%	100%	
8	0%	100%	
9	70%	30%	

2D-LC

Mi n	Sol A: MeOH:H₂O (10:90 v/v) (0.1% FA & 20 mM NH₄HCO₂)	Sol B: ACN:MeOH:IPA (20:30:50 v/v) (0.1% FA & 20 mM NH ₄ HCO ₂)
0	60%	40%
2	40%	60%
4	20%	80%
5	15%	85%
6	10%	90%
7	5%	95%
8	0%	100%
12	0%	100%
13	60%	40%



Figure 1. Due to poor absorbance of the lipid species with the UV detector, total ion chromatograms from MS were used to guide multiple heart cutting 2D-LC



experiments. High resolution experiments with 10 - 6 second cuts were performed across 5 different 1minute regions of the first dimension chromatogram. This resulted in 50 second dimension chromatograms each with a 14 minute gradient which is effectively a comprehensive experiment, but with longer 2D runs than with the typical comprehensive workflow.

2



Figure 3. 2D Extracted chromatograms for selected lipids

IM-MS

lon mobility was also evaluated for lipid separations. Plotting m/z vs. CCS creates a conformational space for the lipid standards. Lipids generally group according to class. Ceramides occupy a larger conformational space than the lyso lipids with similar m/z. The conformational space for the other lipids follows the trend of PC > PE > PG > PS which mimics their head groups in size. Additionally we see lipids with longer tails having larger CCS values. Degree of saturation also plays a role in the trends observed.



Figure 4. Conformational space plot of the lipid standards.

Results and Discussion





Figure 6. Identification results for internal standards spiked in the bovine heart extract are shown in the bar chart. Identifications were based on a database search (DB) and library MS/MS spectral match (Lib) for both positive and negative mode. The pie charts summarize lipid class results for those lipids identified with both DB and Lib. match in both positive and negative mode.

References

¹Unraveling the Complexity of Lipidomes by Multiple Heart Cutting Q-TOF LC-MS with the Agilent 1290 Infinity 2D-LC Solution, G. Vanhoenacker, R. t'Kindt, F. David, P. Sandra, and K. Sandra, Application Note: Biotherapeutics and Biosimilars, 2015.

²An Interlaboratory Evaluation of Drift Tube Ion Mobility Collision Cross Section Measurements, S. M. Stow, T. J. Causon, X. Zheng, R. T. Kurulugama, T. Mairinger, J. C. May, E. E. Rennie, E. Baker, R. Smith, J. A. McLean, S. Hann and J. C. Fjeldsted, Anal. Chem. 2017, 89(17), 9048-9055.

For Research Use Only. Not for use in diagnostic procedures.



This information is subject to change without notice.

© Agilent Technologies, Inc. 2018