

Ultraviolet photodissociation (UVPD) mass spectrometry for structural characterization of lipids in biological matrices on chromatographic time scales

Rahul Deshpande¹, Bashar Amer¹, Susan Bird¹, and Thomas Moehring²

¹Thermo Fisher Scientific, San Jose, USA, ²Thermo Fisher Scientific, Bremen, Germany

ABSTRACT

Purpose

Establish a workflow for structural characterization of lipids using multiple fragmentation techniques including Ultraviolet photodissociation (UVPD).

Methods

Experiments were conducted to characterize lipid structure for resolution of isomers based on sn-position, chain length and double bond positions.

Results

Lipid standard were used for optimization of parameters for characterization of isomers using high resolution chromatographic separation and multiple fragmentation techniques. Resolution of isomers and annotation of unknown lipid species were shown in the bovine liver lipid extract.

INTRODUCTION

Lipidomics is the comprehensive analysis of molecular lipid species. Separation of lipids using liquid chromatography followed by analysis by mass spectrometry has become the method of choice for lipidomic studies. Lipid structures are extremely diverse, in principle rendering over 100,000 unique species by theoretical estimation. The exact number of lipid molecules is even larger due to the presence of structural isomers. The identification of lipids based on their classification is shown in figure 1. Complete structural characterization of complex lipids, such as glycerophospholipids, by mass spectrometry continues to present a major challenge. An ideal analytical tool should be able to characterize the structure of lipids in a discovery platform as well as being feasible for large scale studies. Here we present preliminary work for structural characterization of lipids in a biological matrix on a chromatographic time scale.

FRAGMENTATION

Collision-induced dissociation (CID):

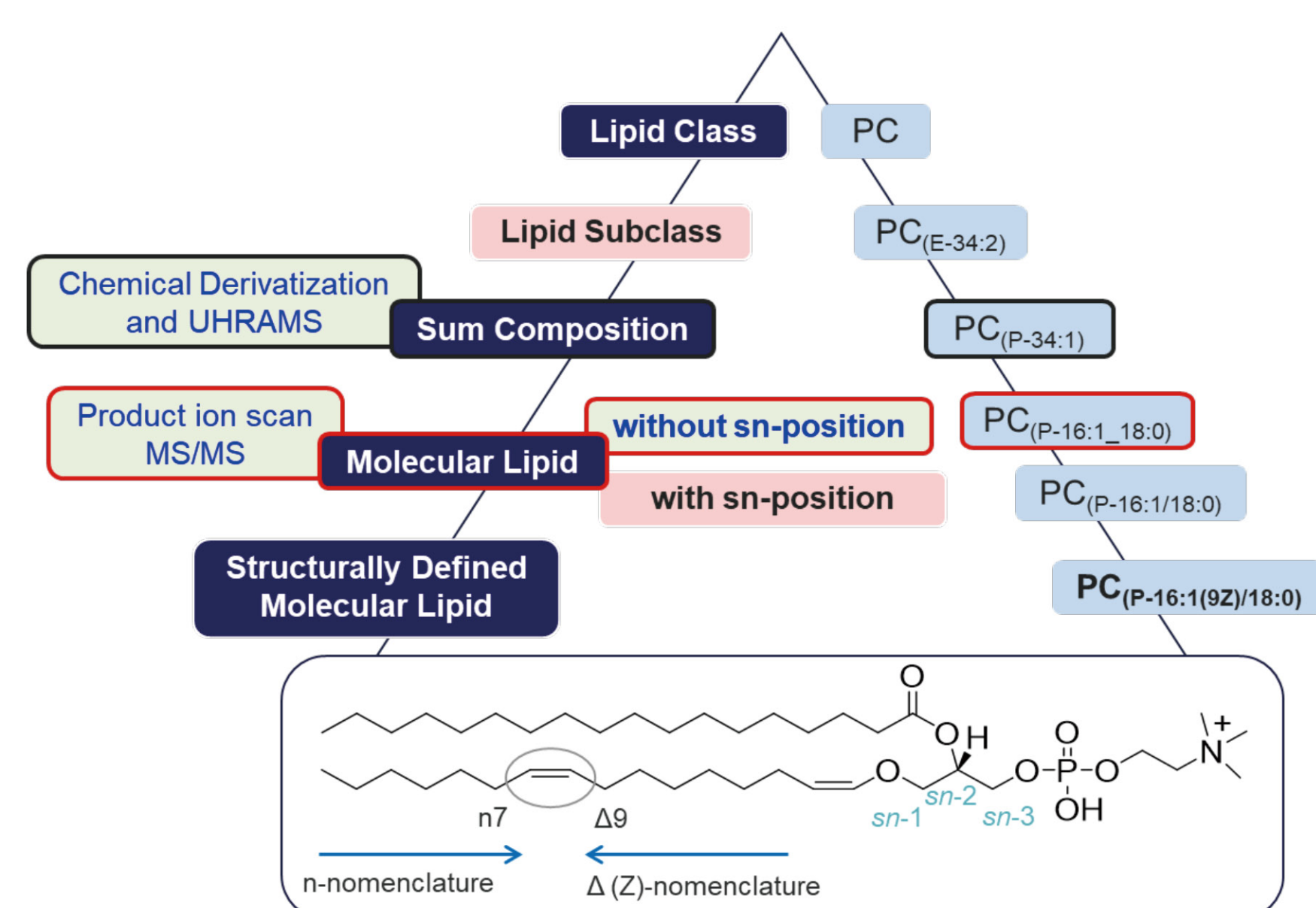
CID or CAD refers to the process whereby the ions kinetic energy has been increased prior to colliding with neutral molecules converting kinetic energy to internal energy that can induce fragmentation. Cannot detect low MW fragments (1/3 Rule)

Higher-energy collisional dissociation (HCD):

HCD is a CID technique associated with Thermo Scientific Orbitrap instruments. Voltage offsets between components increase the kinetic energy of the precursor ion and dissociated in the ion routing multipole cell from collisions with nitrogen molecules. Can detect low mw ions.

Ultraviolet photodissociation (UVPD) -213 nm:

UVPD differs from CID and HCD in that it utilizes photons generated from a laser source to increase the internal energy of a selected precursor ion until there is sufficient internal energy present to overcome the barrier to dissociation. It generates complimentary fragments to HCD/CID.



MATERIALS AND METHODS

Chemicals and Standards

Lipid standards and bovine liver lipid extract were purchased from Avanti Lipids

LC Methods

Column: Thermo Scientific™ Accucore™ C30, 2.6 μm, 150 x 2.1 mm (RP)
Column Temperature: 45 °C
Autosampler Tray Temperature: 15 °C
Injection Volume: 2 μL
Flow rate: 0.26 ml / min
Run Time: 45 Minutes

Mass Spectrometer

Thermo Scientific™ Orbitrap IQ-X™ Tribrid Mass Spectrometer was used.

MS Conditions

MS¹ at 240K resolution (FWHM @ m/z 200) and data-dependent MSⁿ experiments were performed.

Data Processing

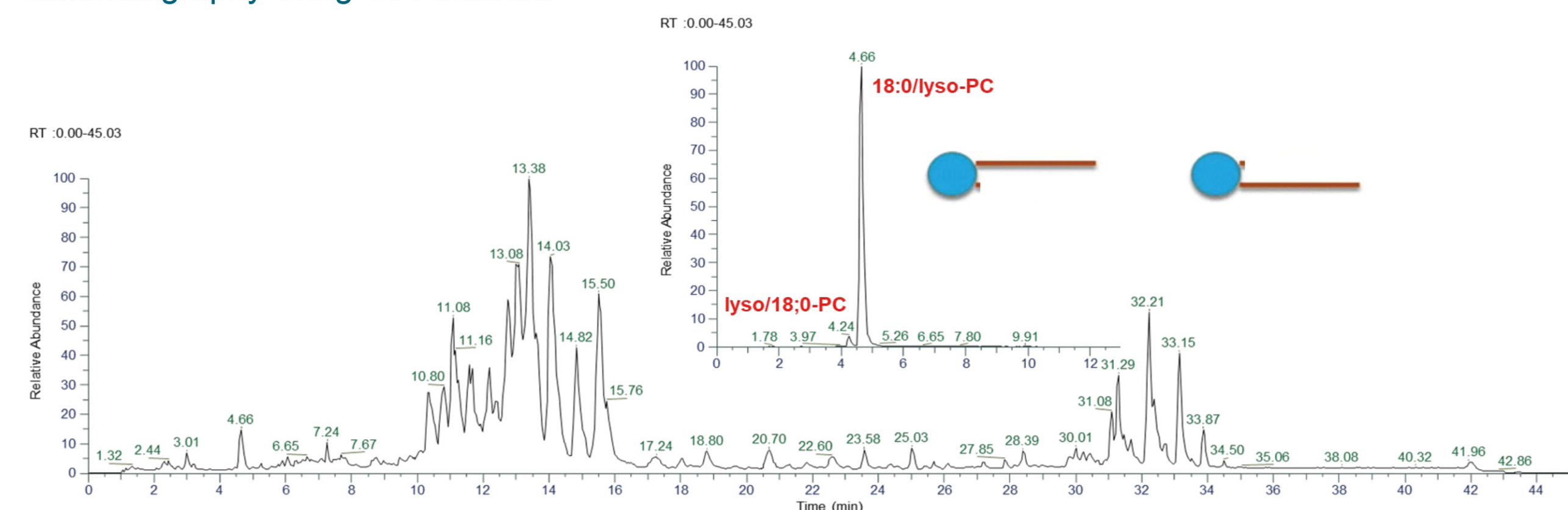
Thermo Scientific™ LipidSearch™ 5 was used for data analysis.

RESULTS

REGIOISOMERS (SN-POSITION)

Lysophosphatidylcholines (lysoPCs) are products or metabolites of phosphatidylcholines (PCs), which are structural components of animal cell membranes. LysoPCs occur in two forms, with the fatty acyl groups at positions 1 (sn-1) or 2 (sn-2) on the glycerol backbone. These two forms can be resolved with high resolution chromatography.

Figure 2 Separation of isomeric lysoPC's in a biological matrix using high resolution liquid chromatography using C30 column.



Phosphatidylcholines (PCs) contain two fatty acyl chains which can result in two regioisomers. These can be resolved using UVPD for MS³ fragmentation of specific neutral loss fragments.

Figure 3 Mass Spectrometric Template on IQ-X Utilizing Complimentary Multiple Fragmentation Strategies for determination and resolution of structural isomers.

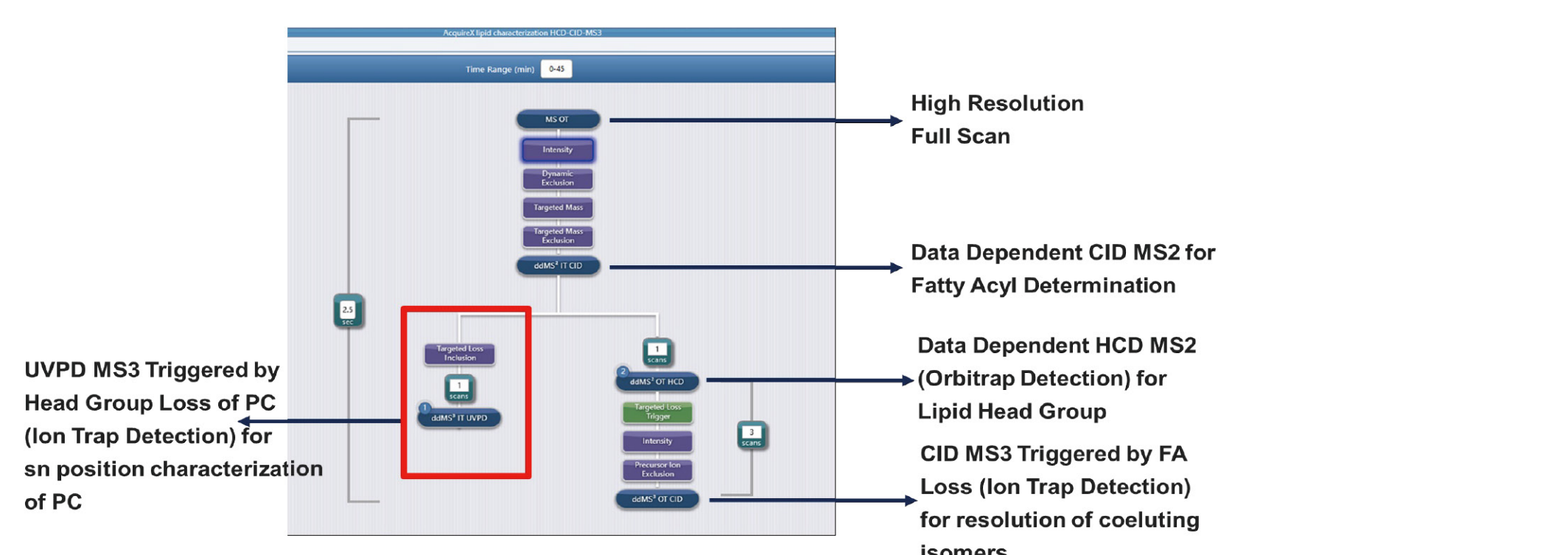
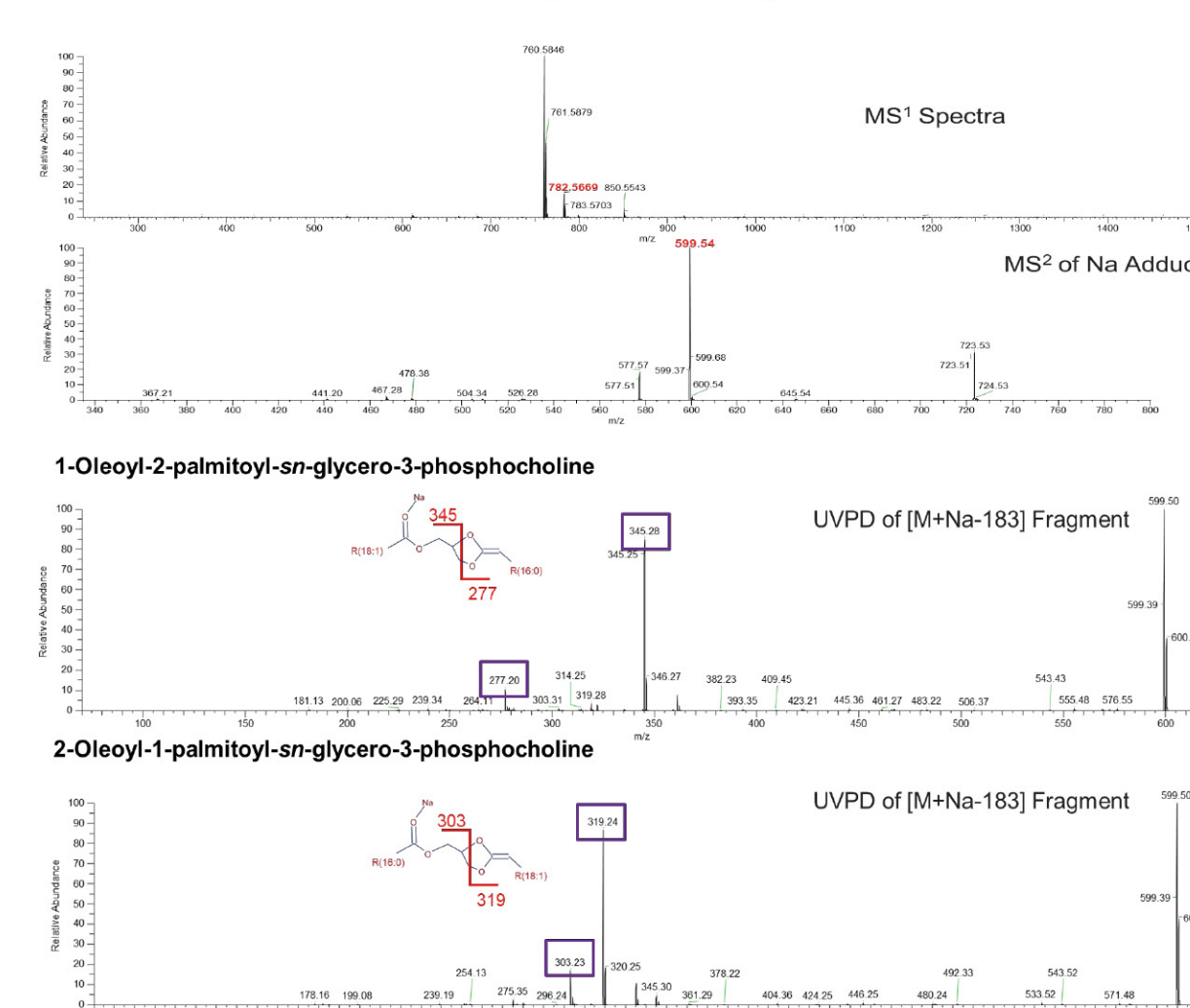


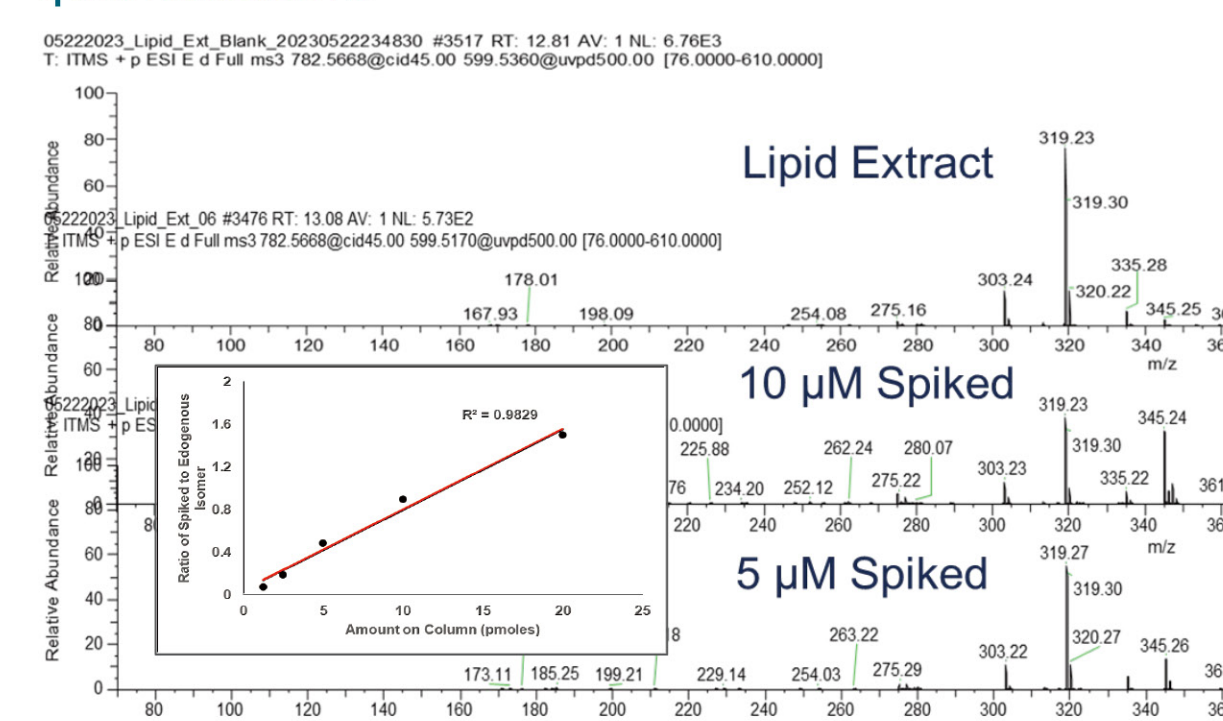
Figure 4 UVPD based MS³ fragmentation for sn-position determination of PC. Lipid standards PC(16:0/18:1) and PC(18:1/16:0) were used for determination of diagnostic fragments for differentiation in biological sample.



QUANTITATION

Bovine Liver Lipid Extract was spiked with 1-Oleoyl-2-palmitoyl-sn-glycero-3-phosphocholine at different ratio. Relative quantification of each isomer was calculated by ratio of diagnostic fragments.

Figure 6 Relative Quantification of regioisomers in biological matrix. Ratio of the diagnostic fragments 345.24 and 319.24 were used for the quantification.



Double Bond

Lipid Standards 1,2-dioleoyl-sn-glycero-3-phosphatidylcholine {PC(18:1(9Z)/18:1(9Z))} and 2-dipetroselinoyl-sn-glycero-3-phosphocholine {PC(18:1(Δ6)/18:1(Δ6))} which are PC isomers differing in the position of the double bond were baseline separated by the high-resolution chromatographic method using C30 column. However, none of these two PC's were present in the bovine liver lipid extract with an isomer at different retention time.

Figure 8 Isomers of PC differing in their double bond position were separated by the high-resolution chromatographic method using C30 column.

