

Differentiating sugar phosphate isomers using high resolution mass spectrometry coupled with MSⁿ, collision induced dissociation, and ultraviolet photodissociation

Brandon Bills¹, Sunandini Yedla¹, Rahul Deshpande¹, Bashar Amer¹, Susan Bird¹, Vlad Zabrouskov¹

¹Thermo Fisher Scientific, 355 River Oaks Pkwy, San Jose, CA, U.S.A., 95134

Abstract

Purpose: Differentiating isomers poses an analytical challenge, especially if they cannot be resolved chromatographically. While tandem mass spectrometry can distinguish structurally distinct isomers based on unique fragments, there isn't always a Higher-energy Collisional Dissociation (HCD) fragment that is diagnostic when the structures are similar. Here we demonstrate how using MSⁿ fragmentation, Collision Induced Dissociation (CID), and Ultraviolet Photodissociation (UVPD) can be used to obtain diagnostic fragments for structurally similar isomers.

Methods: A set of six sugar phosphate isomers: glucose-1-phosphate, galactose-1-phosphate, glucose-6-phosphate, galactose-6-phosphate, mannose-6-phosphate and fructose-6-phosphate were analyzed individually by LC-MS. Spectra were collected using HCD, CID, MSⁿ, and UVPD then evaluated for diagnostic fragments.

Results: Each of the sugar-phosphate standards exhibited at least one fragment unique to that isomer at that RT.

Introduction

Sugar phosphates, sugar molecules with a covalently bound phosphate, are an important intermediate in carbohydrate metabolism and a building block for oligonucleotides. This class of compounds can be structurally very similar with poorly resolved chromatograms and similar mass spectra making characterizing what compounds are present difficult. In this work we use a range of different fragmentation options available on the Thermo Scientific™ Orbitrap™ IQ-X™ Tribrid™ mass spectrometer to investigate individual sugar phosphate standards to identify unique fragments that are diagnostic of a specific compound.

Materials and methods

Sample Preparation

Each sugar (glucose-1-phosphate, glucose-6-phosphate, fructose-6-phosphate, mannose-6-phosphate, galactose-1-phosphate and galactose-6-phosphate from Sigma-Aldrich) was prepared at 1 mM in 60:40 Acetonitrile:Water as well as a mixture of all six sugar standards at 0.5 mM in quenched plasma (plasma after an acetonitrile protein crash).

Test Method

LC-MS: All samples were run on a Thermo Scientific™ Vanquish Horizon™ LC system using a ZIC - pHILIC column (Sigma Aldrich) with a mobile phase of A : 5mM Ammonium carbonate+0.1% NH₄OH in 100% Water, B : 100% Acetonitrile using a Gradient of 80% ACN to 20% in 20 mins. The detector was a Thermo Scientific Orbitrap IQ-X Tribrid mass spectrometer run in positive ion mode using DDA MSⁿ, HCD, CID, and UVPD fragmentation.

Data Analysis

Chromatograms and mass spectra were processed using Thermo Scientific™ Freestyle™ software. Structural predictions of the fragments were generated using Thermo Scientific™ Mass Frontier™ software.

Figure 1. Thermo Scientific Horizon Vanquish LC and Orbitrap IQ-X Tribrid mass spectrometer.

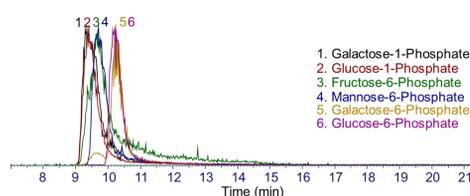


Results

Chromatography

Running each standard individually, it was found that four of the six sugar phosphates eluted with significant chromatographic overlap while the remaining two eluted at a later retention time. In each case, a peak at for the protonated (m/z 261.0370) and sodiated adducts (m/z 283.0189) was observed.

Figure 2. Overlaid XIC chromatogram for the six protonated sugar phosphate standards.



Ultraviolet Photodissociation diagnostic peaks

UVPD spectra for each standard were investigated for peaks that were diagnostic, present for only one standard at that retention time. The first four standards were compared for similar peaks as were the last two that eluted after 10 minutes.

Figure 3. Diagnostic UVPD fragment for protonated adduct of galactose-1-phosphate.

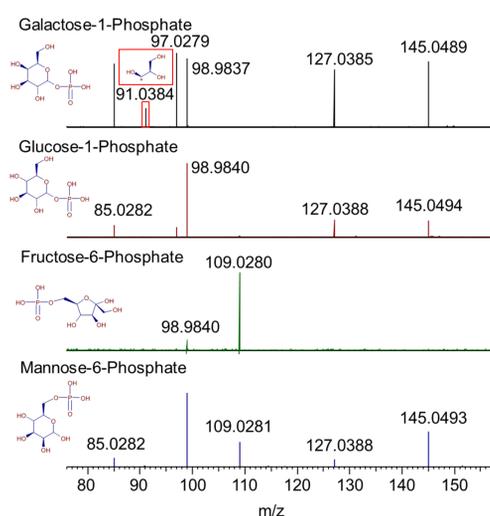


Figure 4. Diagnostic UVPD fragment for sodiated adduct of glucose-1-phosphate. The relevant section of the spectra has been scaled up for visibility.

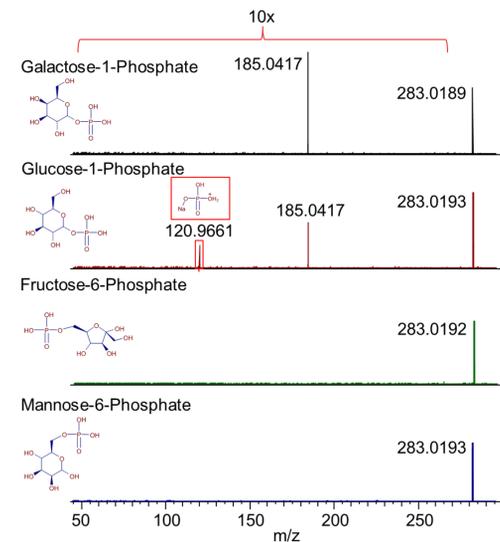
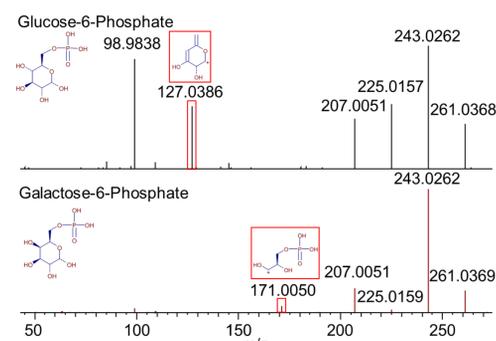


Figure 5. Diagnostic UVPD fragments for proton adduct of glucose-6-phosphate and galactose-6-phosphate.



MSⁿ CID diagnostic peaks

HCD MS² and CID MS³ spectra for each standard were investigated for peaks that were diagnostic for fructose-6-phosphate and mannose-6-phosphate

Figure 6. Diagnostic HCD MS² fragments for sodium adduct of fructose-6-phosphate and mannose-6-phosphate.

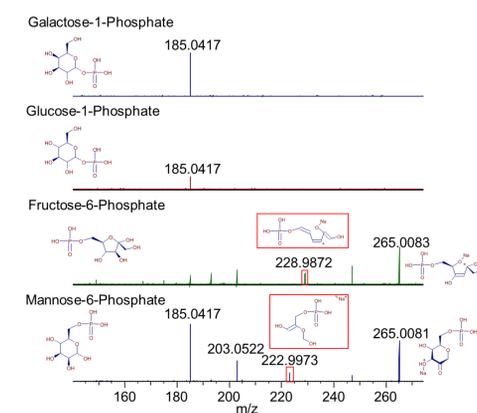


Figure 7. Diagnostic CID MS³ fragments of the MS² precursor at m/z 265.0081 from mannose-6-phosphate.

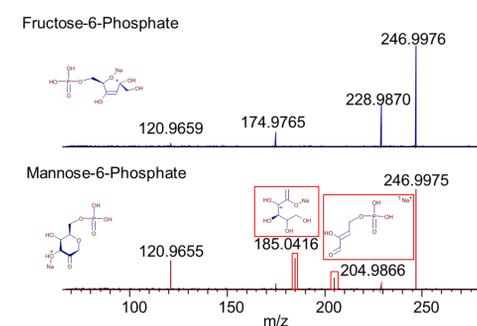


Figure 8. Diagnostic UVPD MS² fragment collected in the ion trap for sodium adduct of glucose-1-phosphate. The region of importance has been scaled up for visibility.

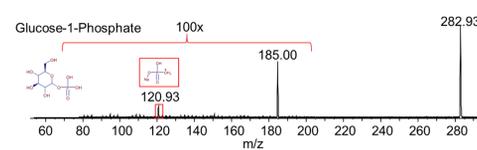
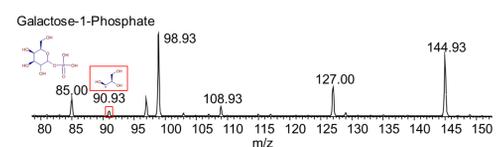


Figure 9. Diagnostic UVPD MS² fragment collected in the ion trap for proton adduct of galactose-1-phosphate.



Conclusions

Six isomeric sugar phosphate compounds were investigated using HCD, MSⁿ, CID, and UVPD fragmentation for diagnostic fragments

- Four coeluting isomers as well as a second set of two coeluting isomers could be clearly differentiated using the fragmentation options available on the Orbitrap IQ-X Tribrid mass spectrometer
- Even when mixed together in a complex matrix, these diagnostic fragments were visible using the high sensitivity of the ion trap

Acknowledgements

We would like to thank Hardik Shah at the University of Chicago Medicine Comprehensive Cancer Center, Chicago, IL, USA and Rohit Sharma at the Department of Molecular Biology, Massachusetts General Hospital, Harvard Medical School, Boston, MA, USA.

Trademarks/licensing

© 2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

PN: PO66162-EN0422S