

Creating Unique Fragmentation: Small Molecule Structural Elucidation Using UVPD

Jean-Jacques Dunyach, Scott Peterman, Seema Sharma, Romain Huguet, Vlad Zabrouskov, Christopher Mullen, Thermo Fisher Scientific, San Jose, CA, USA

ABSTRACT

Purpose: Utilizing ultraviolet photodissociation (UVPD) to increase the fragmentation efficiency resulting in more structurally informative product ions relative to high energy collisional dissociation (HCD).

Methods: Perform comparative LC-MS/MS experiments on a series of standards as well as a set of hepatocyte metabolite incubations using 213 nm laser as compared to beam-type fragmentation.

Results: For all standards evaluated, compounds with more than 3 Pi bonds showed greater fragmentation using UVPD than HCD as well as resulting in many more unique fragments with both even and odd electrons. In addition, UVPD provides greater diagnostic fragmentation to localize the site of conjugation by retaining the entire glucuronide modification.

INTRODUCTION

Small molecule structural elucidation using LC-MSⁿ presents significant challenges due to the overwhelming structural diversity, lack of mobile protons and/or multiple basic/anionic sites, and biotransformations. Collisional activation/dissociation-based methods results in preferential unimolecular dissociation pathways that may limit the formation of structurally informative product ions. This issue is particularly common for conjugated metabolites or fused-ring structures. We have developed instrument methods designed to leverage the commercially available 213 nm UVPD source on the Thermo Scientific™ Orbitrap Tribrid™ mass spectrometers to increase the structural elucidation capabilities for a series of small molecule standards and a set of isobaric conjugated metabolites from Phase II drug incubations. The LC-MSⁿ method consisted of standard data dependent acquisition (DDA) and dynamic exclusion (DE) settings changing only the dissociation methods between HCD and UVPD. Evaluation of the number of unique product ions and structural coverage were used to determine the effectiveness of UVPD relative to collisional activation, as well as gauge the limitations of UVPD based on the compound structure.

MATERIALS AND METHODS

Sample Preparation

Two sets of standards were used to initially evaluate UVPD performance and optimal settings. One standard contained an equal molar mixture of a range of compounds with various numbers of aryl rings and double bonds. The second set of standards consisted of base structures and its isomeric metabolites such as Olmesartan and isomeric glucuronides structures. Each compound individually analyzed by HCD and UVPD with a common LC gradient and then mixed in an equal molar mixture. Lastly, a set of Phase II hepatocyte incubations of verapamil was analyzed by the comparative methods.

Test Method(s)

All samples were evaluated on a Thermo Scientific™ Orbitrap Fusion™ Lumos™ Tribrid™ mass spectrometer and Thermo Scientific™ Orbitrap Eclipse™ Tribrid™ mass spectrometer using a common LC gradient and DDA/DE method. A binary solvent system was used comprised of A) 0.1% formic acid and B) MeCN with 0.1% formic acid and an 8-minute gradient (10-35%) on a Hypersil Gold aQ column with dimensions of 100 x 2.1 and 1.9 μm particles. The DDA/DE method consisted of a full scan MS (m/z 200-800) and a top 5 method with the MS and MS/MS spectra acquired with a resolving power of 60k and 30k, respectively. The maximum ion fill time for all MS/MS spectra was 54 msec. For HCD activation a normalized collision energy (NCE) of 35% was used while the UVPD pulse duration was maintained at 75 msec.

Data Analysis

All data was processed using Thermo Scientific™ FreeStyle™ Software and Thermo Scientific™ Mass Frontier™ 8.0 spectral interpretation software. Chemical structures were created and saved as a library file as well as used predict product ions using the following settings in Mass Frontier 8.0 software: general fragmentation libraries, both even and odd electron considerations, and a maximum of 5 steps for a unimolecular reaction coordinate. A mass measurement error of 2 ppm was used for initial screening to determine product ions.

RESULTS

Leveraging the unique capabilities of the Tribrid mass spectrometer

The goal of the study was to introduce alternative fragmentation mechanisms to generate complimentary unimolecular dissociation pathways relative to the standard beam-type activation (HCD) to enhance small molecule structural characterization and elucidation workflows. The primary method used in this study was UVPD due to the energy imparted per absorbed photon. A representative energy diagram is presented in Figure 1 comparing common activation methods used on commercial mass spectrometers.¹ Clearly the utilization of 213 nm photons imparts the greatest increase of internal energy per unit of time resulting in altering fragmentation methods from primarily kinetically controlled dissociation pathways to thermodynamically controlled pathways resulting in a higher probability of generating unique and structurally definitive product ions being measured. The Tribrid instrument architecture is ideal for implementing UVPD as the ion flight path and optics enables optimal precursor isolation (quadrupole mas filter), collection (ion routing multipole), trapping for irradiation and collection of the resulting unreacted precursors and all product ions (linear ion trap), and high resolution accurate mass (HRAM) measurement of all ions (Ultra-high field Orbitrap mass analyzer) is controlled by on-the-fly decisions (Dynamic Scan Management) enabling LC-MS/MS analysis on UHPLC time scale (6 second wide peaks) for both high and low-abundant precursor ions. The UVPD laser is contained within the footprint of the instrument to minimize required laboratory space and controlled within the standard instrument control software.

Figure 1. Cartoon representation of relative energy deposition resulting from comparative ion activation methods. The reaction coordinate demonstrates the required energy needed to access various unimolecular dissociation channels resulting in various fragments demonstrating the significant increase in energy deposition following the absorption of a single UV photon. Pi bonds (C=C, C=N) have strong absorption bands overlapping with the commercially available UVPD laser to promote fragmentation.

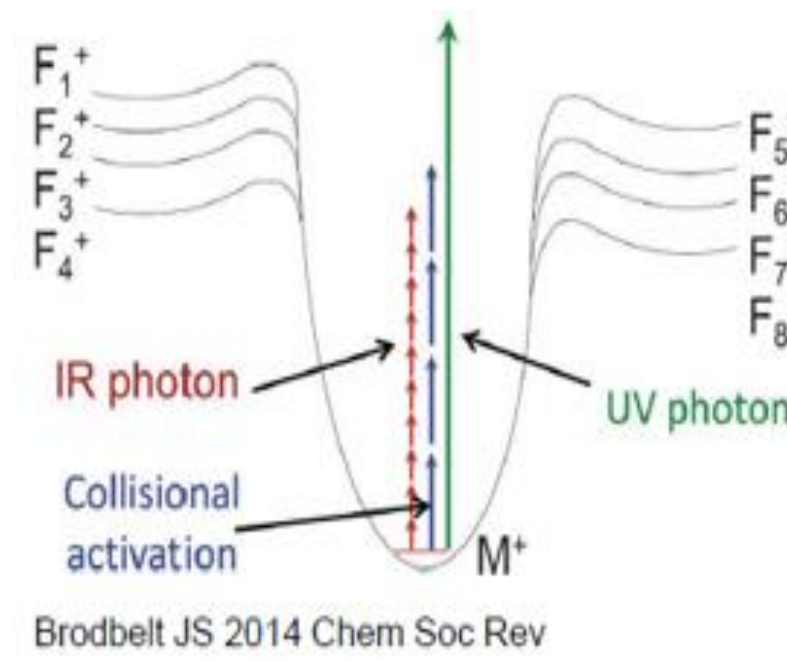


Figure 2. Schematic representation of the Orbitrap Tribrid mass spectrometer with UVPD laser location and introduction into the instrument through the linear ion trap.

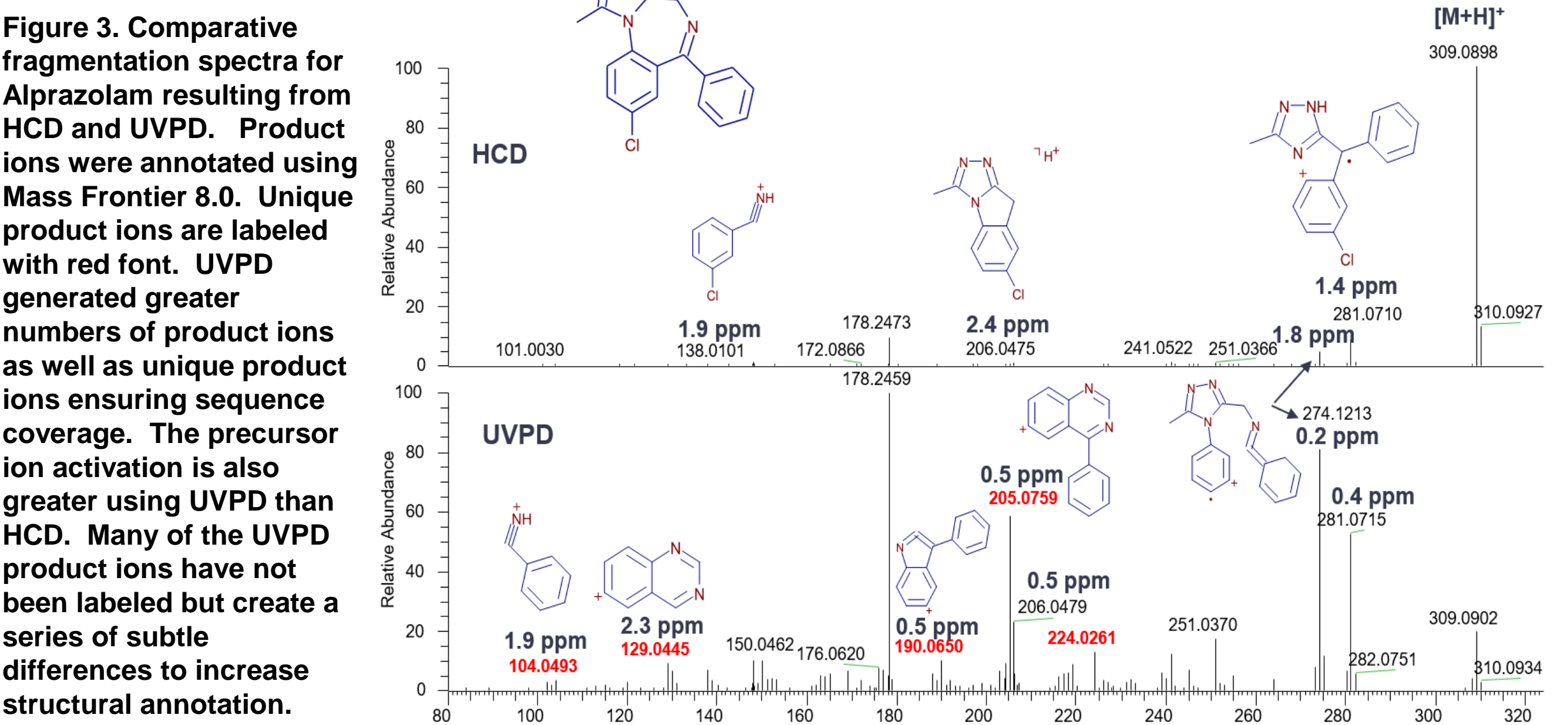
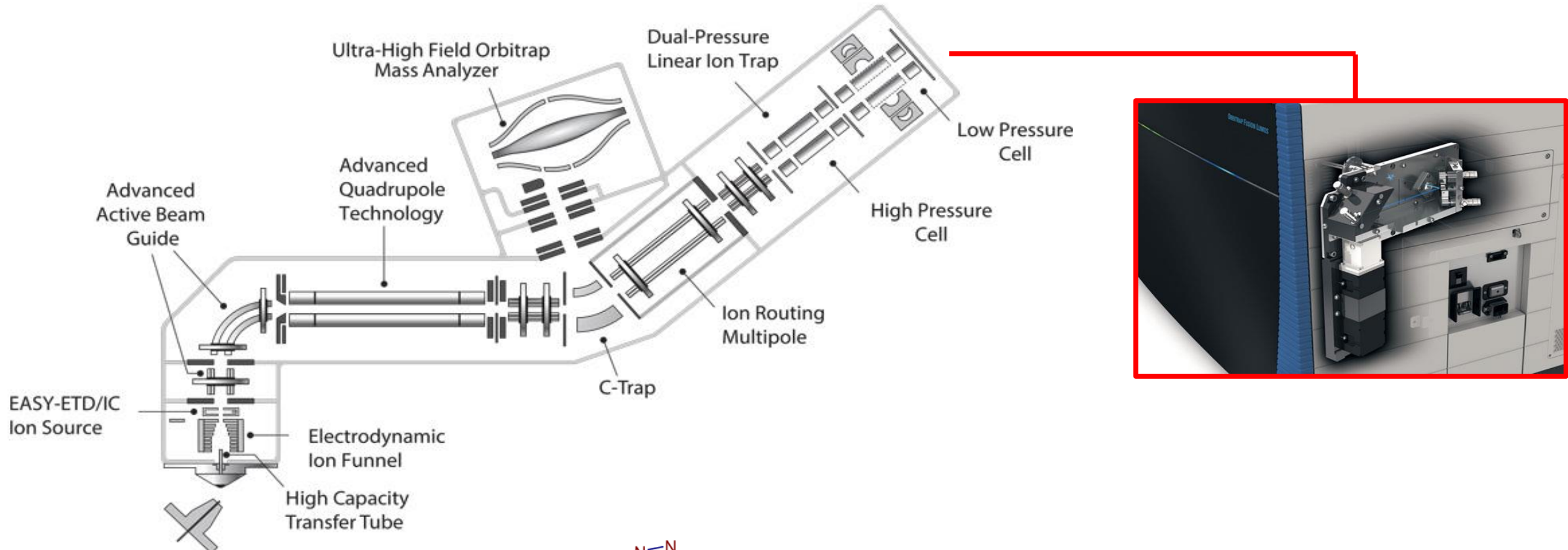


Figure 4. Evaluation of structural elucidation for hydroxy-diclofenac structural isomers using (Fig. 4A) HCD and (Fig. 4B) UVPD. The two isomers shown below demonstrate the challenges using collisional activation of small molecules containing one primary basic site resulting in primarily two fragments that do not enable differentiation. Figure 4B shows the resulting fragmentation following UVPD with much greater fragmentation, and as highlighted, many more unique fragments that are unique per structure that are highlighted in red. Specifically, the highlighted product ion mass range is shown to the right, further demonstrating the importance of high resolution and mass accuracy for detection of key product ions despite it being formed in the 1+ charge state.

