# Implementation of 213 nm Ultra Violet Photodissociation (UVPD) on a Modified Orbitrap Fusion Lumos

Christopher Mullen<sup>1</sup>; Chad R. Weisbrod<sup>2</sup>; Eugene Zhuk<sup>1</sup>; Romain Huguet<sup>1</sup>; Jae C. Schwartz<sup>1, 1</sup>Thermo Fisher Scientific, 355 River Oaks Parkway, San Jose, CA 95134, <sup>2</sup>National High Magnetic Field Laboratory, 1800 E. Paul Dirac Dr., Tallahassee, FL 32310

### **OVERVIEW**

Details of the implementation of UVPD using a 213 nm laser into the Thermo Scientific™ Orbitrap Fusion™ Lumos™ platform are presented, along with results pertinent to top down proteomic experiments.

#### INTRODUCTION

UVPD has been demonstrated to be a useful tool for the analysis of peptides and proteins. Numerous studies have concluded that UVPD promotes all three types of backbone cleavages (resulting in a,b,c,x,y, and z type ions), and that UVPD preserves labile PTMs during activation such as phosphorylation, sulfation, and glycosylation.¹ While its application to the study of large polypeptides and proteins is relatively new, interest in the technique is rapidly increasing due to the sequence coverage provided, advances in proteome fractionation, and top-down informatics.².³ Last year Weisbrod et. al. presented the performance characteristics of UVPD using the 5<sup>th</sup> harmonic of a Nd:YAG laser.⁴ This presentation is a continuation of that initial work and describes in detail the implementation of the 213 nm laser on a modified Orbitrap Fusion Lumos instrument including the performance related to top down protein sequencing.

## **MATERIALS AND METHODS**

#### Sample Preparation

Apomyoglobin from equine skeletal muscle, lyophilized powder, Sigma-Aldrich A8673. 1 pmol/µL in 50/50 MeOH/H<sub>2</sub>0 w/ 0.1% Formic Acid.

Carbonic Anhydrase II from bovine erythrocytes, lyophilized powder, Sigma-Aldrich C2522. 1 pmol/µL in 50/50 MeOH/H<sub>2</sub>0 w/ 0.1% Formic Acid.

#### Test Method(s)

The sequence coverage reported for apomyoglobin and carbonic anhydrase was attained by infusing a 1 pmol/uL solution of the analyte at 3 uL/min into the instrument. The scan parameters were 5e5 precursor agc target, 120K OT resolution, 100 uScans averaged, with the instrument in the low pressure mode (IRM = 3.0 mTorr). The UVPD irradiation time, and hence the number of laser pulses to which the analyte was exposed was varied to find the maximum in sequence coverage.

#### Data Analysis

MS/MS spectra were deconvoluted using Hardklör⁵, and searched against their respective sequence using ProSight Lite<sup>6</sup> with a 10 ppm mass tolerance.

Figure 1. Modified Orbitrap Fusion Lumos mass spectrometer with laser assembly interfaced to the rear of the ion trap manifold. Shown is the assembly flange, laser head, steering optics, and vacuum interface window which are all behind the side cover.



## **RESULTS - IMPLEMENTATION**

#### **Laser Description**

The laser employed is a Q3 series passively Q-switched Nd:YAG laser (CryLaS GmbH) outputting the 5<sup>th</sup> harmonic at 213 nm. The laser has an average power consumption of approximately 20 Watts, and is passively cooled on a heat sink that also functions as its mounting bracket. The laser pulse characteristics and relative size can be seen in Table 1 and Figure 2, respectively.

 Table 1. Laser Characteristics

	Reference Value	Unit
Output Power (quasi cw)	3.75 ±0.5	mW
Pulse Energy	$1.5 \pm 0.2$	μJ
Peak Power	1.5	kW
Pulse Rep Rate	2.5	kHz
Pulse Width	< 1	nsec
Beam Dia	450 ±200	μm
•		

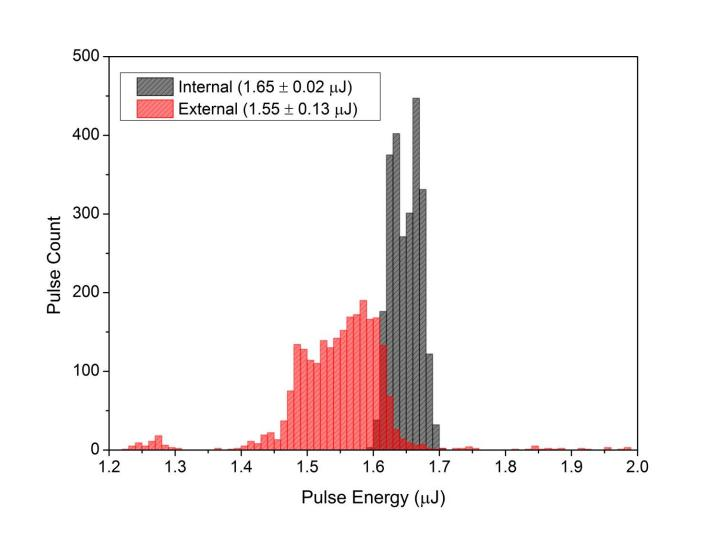
Figure 2. CryLaS 213 nm Nd:YAG laser



The laser pulse energy is  $1.5 \pm 0.2$  uJ/pulse at 2.5 kHz repetition rate (0.4 msec/pulse). The beam diameter including divergence at the center of the ion trap is slightly larger than the simulated ion cloud diameter at normal agc targets, and therefore no focusing optics are required. The laser trigger is provided by the instrument as is synched to the scan function. Photoactivation occurs in the low pressure cell of the dual cell quadrupolar linear ion trap, while m/z analysis can occur in either the linear trap or the Orbitrap mass analyzer.

The laser's pulse to pulse stability is very good and has been optimized for variable delays between pulse trains to account for the overhead between scans associated with manipulating ions and m/z analysis. Figure 3 shows the pulse variability for 2500 pulses measured using the CryLaS internal trigger and with the external trigger supplied by the instrument during the scan function.

Figure 3. Histogram of the laser pulse energy measured using a Coherent J-10MT-10kHz-USB power meter. Black trace is from internal triggering of the laser at 2.5 kHz. Red trace is external triggering of the laser using the instrument scan function with 125 pulses/50 ms activation time and an approximately 50 msec pause between pulse trains.



#### **UVPD Laser Integration to the Orbitrap Fusion Lumos mass spectrometer**

As can be seen in Figure 4, the 213 nm laser is affixed to the system via a modified ion trap chamber back flange, which also houses the steering optics and vacuum interface window. All optics within the laser path are high quality UV grade fused silica, and are appropriately coated.

Figure 4. UVPD module and interface to the dual cell linear ion trap. [1] – Laser heat sink, [2] – Mirror1, [3] – Mirror2, [4] – Vacuum flange, [5] – fused silica vacuum interface window, [6] – dual cell linear ion trap assembly (sectioned).

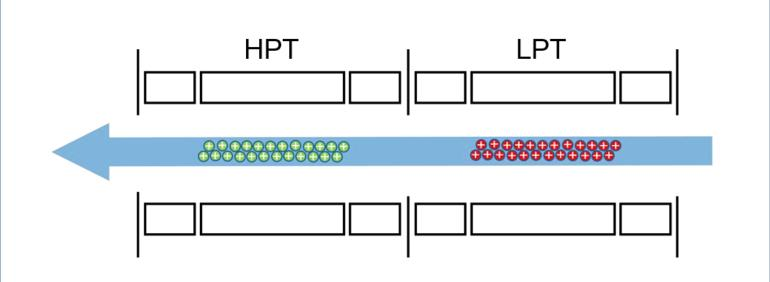


The laser beam is manually aligned to the ion cloud in both ion traps using the actuators at each of the two mirrors. Alignment to the ion clouds in both ion traps was deemed advantageous as it ensures the maximum overlap of the laser beam with the ion cloud while minimizing the chance for photo-generated noise in the system. The Alignment Quality Score (AQS) characterizes the quality of the overlap and is described as:

$$AQS = \frac{\%PD_{HPT}}{\%PD_{HPT} + \%PD_{LPT}}$$

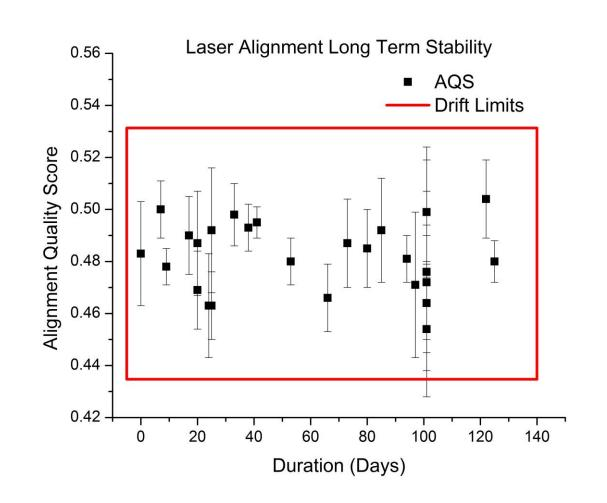
where %PD is the % photodissociation observed for a fixed amount of activation time, and HPT and LPT are the high pressure and low pressure cell, respectively. We therefore strive for an AQS = 0.5, while simultaneously maximizing the %PD in each cell, Figure 5.

Figure 5. Depiction of good laser alignment. The laser beam interacts with the total ion population in both ion traps.



The UVPD module, once aligned, shows good stability with time, leading to a robust implementation requiring little additional user intervention, Figure 6.

Figure 6. Laser beam – ion cloud overlap long term stability.



## RESULTS – TOP DOWN PERFORMANCE

#### **Sequence Coverage**

The sequence coverage for UVPD of apomyoglobin and for carbonic anhydrase as a function of activation time, and for various charge states, is presented in Figure 7. The data in the figure was acquired according to the conditions described in the method section. In general, at the optimum activation times we achieve approximately 55%, and 30% sequence coverage, respectively for 100 uScans averaged. Figure 8 compares the total bond coverage for various dissociation methods, as well as the combined result of 95% which indicates the unique sequence information provided by UVPD. The data in Figure 8 represents a compilation of multiple activation times for each activation method. They are, 5,10, 17.7, 25, and 35 msec for UVPD, 3, 4, and 5 msec for ETD, 35 %NCE for CID and 10 and 12 %NCE for HCD.

## Figure 7. Sequence coverage vs. UVPD activation time for Apomyoglobin (16.8 kDa) and Carbonic anhydrase (29 kDa) obtained using the conditions described in the methods section for various charge states.

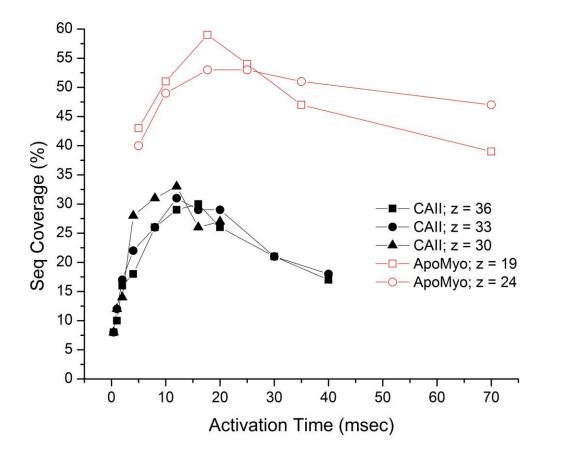
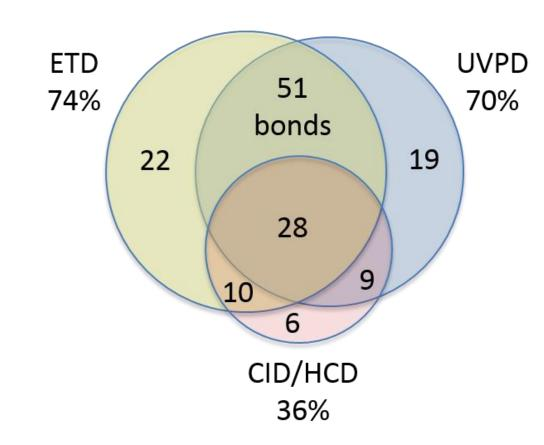


Figure 8. Bond coverage comparison for apomyoglobin [M+24H]<sup>24+</sup> charge state for UVPD, ETD, CID, and HCD, and for the combination.



## **CONCLUSIONS**

- A compact, robust 213 nm solid state laser has been incorporated into the Orbitrap Fusion Lumos platform for the routine implementation of UVPD.
- Details of the implementation, including the laser characteristics, alignment strategy and stability, are described.
- The 213 nm UVPD performance for intact protein standards is shown to be very good, indicating
  its utility for top-down applications, including biotherapeutics. (Please see Fornelli, TP087).

## **REFERENCES**

- 1. Broadbelt JS. "Ion Activation for Peptides and Proteins" Anal Chem 2016;88:30-51.
- Shaw JB, Li W, Holden DD, Zhang Y, et. al. "Complete protein characterization using top-down mass spectrometry and ultraviolet photodissociation" J. Am. Chem. Soc. 2013;135:12646-12651.
- 3. Cleland TP, DeHart CJ, Fellers RT, et.al. "High-throughput analysis of intact human proteins using UVPD and HCD on an Orbitrap mass spectrometer" *J. Proteome Research* 2017;15(5):2072-2079
- 4. Weisbrod CR, Zhuk E, Dunyach JJ, and Schwartz, JC. "Performance considerations for ultraviolet photo-dissociation using the Nd:YAG 5<sup>th</sup> harmonic (213 nm)" ASMS 2016; ThOB 10:10am.
- 5. Hoopman MR, Finney GL, and MacCoss MJ. "High-speed data reduction, feature detection, and MS/MS spectrum quality assessment of shotgun proteomics data sets using high-resolution mass spectrometry" *Anal Chem* 2007:Aug 1;79(15):5620-5632.
- 6. Fellers RT, Greer JB, Early BP, Yu X, LeDuc, RD, Kelleher NL. "ProSight Lite: Graphical software to analyze top-down mass spectrometry data" *Proteomics* 2015: Feb 26;15(7):1235-1238.

#### TRADEMARKS/LICENSING

© 2017 Thermo Fisher Scientific Inc. All rights reserved. ProSight is a trademark of Proteinaceous, Inc. All other 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.