

# Performance Evaluation of the Q Exactive HF Hybrid Quadrupole-Orbitrap Mass Spectrometer for High-Throughput Top-Down Proteomics

Eugen Damoc,<sup>1</sup> Ping Yip,<sup>2</sup> Leena Valmu,<sup>3</sup> Alexander Cherkassky,<sup>2</sup> Bernard Delanghe,<sup>1</sup> Eduard Denisov,<sup>1</sup> Helene Cardasis,<sup>2</sup> Jason Neil,<sup>2</sup> Alexander Makarov,<sup>1</sup> Jim Stephenson<sup>2</sup>  
<sup>1</sup>Thermo Fisher Scientific, Bremen, Germany; <sup>2</sup>Thermo Fisher Scientific, Cambridge, MA, USA; <sup>3</sup>Thermo Fisher Scientific, Vantaa, Finland

## Overview

**Purpose:** Evaluation of the Thermo Scientific™ Q Exactive™ HF hybrid quadrupole-Orbitrap mass spectrometer for high-throughput top-down proteomics.

**Methods:** Top-down analysis of an *Escherichia coli* extract using the data dependent “TopN” method with and without chromatographic separation.

**Results:** We demonstrate utility and applicability of the Q Exactive HF mass spectrometer to perform high-throughput top-down proteome analysis.

## Introduction

Major goals in every top-down proteomics experiment are protein identification and characterization. The strategy used to achieve these goals involves high-resolution mass measurement of intact protein ions followed by their fragmentation and analysis in the mass spectrometer. In spite of enormous improvements in terms of speed and sensitivity in FTMS instrumentation over the last few years, top-down LC-MS/MS in large scale proteome analyses will further benefit if high resolution analysis at higher detection speed would be possible. Furthermore, improvement to the current generation of charge assignment and protein deconvolution algorithms to handle complex top-down data will lead to more efficient, complete, and accurate protein identification. Here we demonstrate the improved performance of the Q Exactive HF hybrid quadrupole-Orbitrap mass spectrometer in a series of high-throughput top-down proteomics experiments in conjunction with a new algorithm for charge assignment and protein deconvolution. Furthermore, a multiplex SIM approach to isotopically resolve multiple charge states of proteins up to 50 kDa at LC timescale is presented.

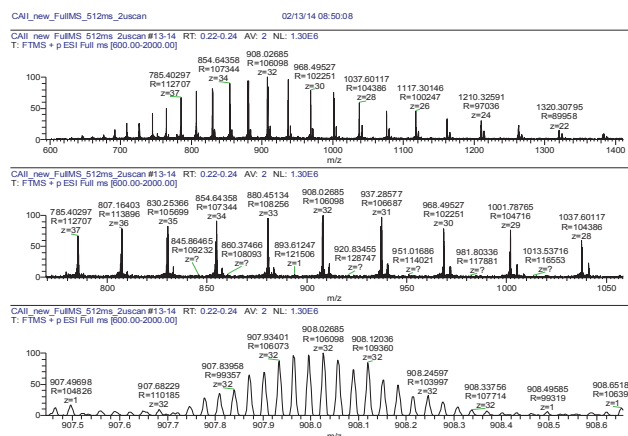
## Methods

Direct infusion experiments using intact carbonic anhydrase II were carried out to evaluate the ability of the Q Exactive HF instrument to perform top-down analysis. Also, top-down microbial proteome analysis was performed by LC-MS/MS or direct static nanospray utilizing an *E. coli* extract. 1–2 µg of protein sample was loaded onto a Thermo Scientific™ PepSwift™ Monolithic PS-DVB (200 µm × 25 cm) EASY-Spray™ column, and four different LC gradients (5, 15, 30, and 60 min) were run on a Thermo Scientific™ EASY-nLC™ 1000 system. A data-dependent “Top-N” method using the “high-high” approach was employed to deliver high resolution and high mass accuracy in both MS and MS/MS modes, using the Q Exactive HF hybrid quadrupole-Orbitrap mass spectrometer. Proteoforms were identified using a new charge assignment and protein deconvolution algorithm. Furthermore, the high-throughput top-down proteomics data was analyzed using Thermo Scientific™ ProSightPC 3.0 software. Multiplexed SIM experiments were performed using LC/MS analysis of intact enolase.

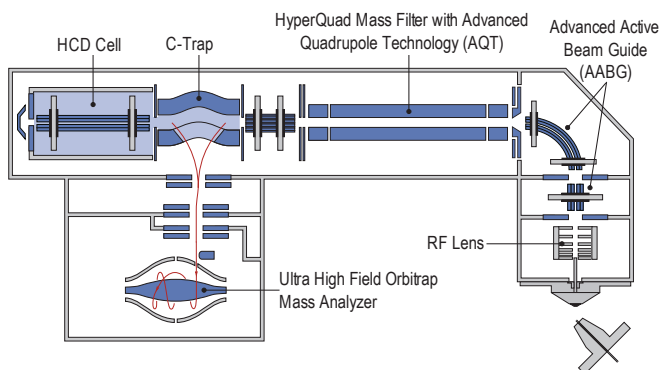
## Results

With the implementation of the compact ultra-high field Thermo Scientific™ Orbitrap™ analyzer on the Q Exactive HF instrument (see Figure 1), the resolving power has been increased by 1.8 fold over that of the previous Orbitrap detector. This enables high-resolution analysis at high detection speed which makes the HF instrument more suitable for top-down analysis at LC time scale. The novel Intact Protein Mode allows adjustment of the trapping gas pressure and optimizes the control logic of the instrument to analyze intact proteins with masses up to 50 kDa with isotopic resolution. Carbonic anhydrase II with a molecular mass of 29 kDa was used to evaluate the ability of the Q Exactive HF instrument to perform top-down analysis. Figure 2 shows results of an experiment, where full MS scans were recorded at a resolving power setting of 240,000 (FWHM at  $m/z$  200) and AGC target value of 3e6. The figure shows an averaged spectrum over 2 seconds, where the isotopes are baseline resolved and the charge states are properly assigned. Figure 3 is retrieved from a second experiment, where an MS/MS scan with higher-energy collisional dissociation (HCD) of the charge state 34+ at a collision energy of 20 eV was performed. The AGC target value was 1e6 at a resolving power setting of 120,000 (FWHM at  $m/z$  200) with 4 µscans in 1 second acquisition time. 36 *b*-type and 28 *y*-type fragment ions were identified using ProSight PC 3.0 software.

**FIGURE 2. Full-MS spectrum of intact carbonic anhydrase II (2 × 2 µscans @ 240k res. pwr. → acq. time: 2 seconds) with baseline resolution of the isotopic pattern.**



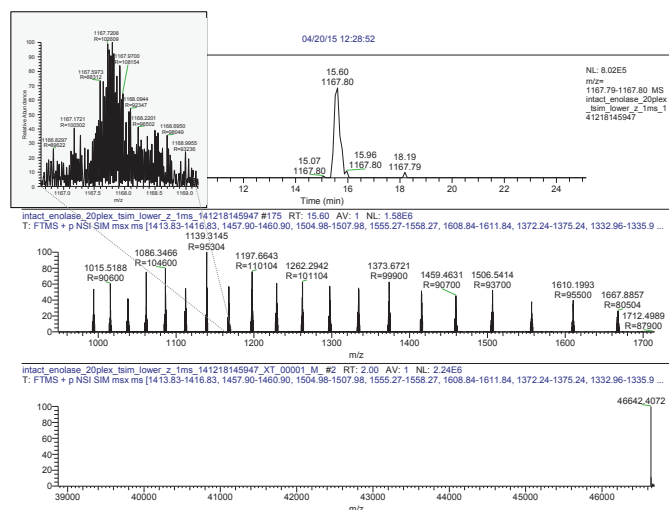
**FIGURE 1. The Q Exactive HF instrument layout.**





Most of the proteins identified using "Top-N" "high-high" method have molecular weights < 35 kDa, which is why a multiplex SIM approach was tested to see whether the mass range of isotopically resolved proteins can be extended beyond this limit. With this approach, different charge states of the same protein or of different proteins can be first selected using the quadrupole, then trapped in the HCD cell, and detected all together with the Orbitrap analyzer. Using intact enolase we could demonstrate that proteins up to about 50 kDa can be analyzed with isotopic resolution at LC time scale (see Figure 7).

**FIGURE 7. Multiplex SIM spectrum of 20 consecutive charge states of intact enolase (10 μscans @ 240k res. pwr. → acq. time: ~ 5 seconds).**



## Conclusion

- Q Exactive HF mass spectrometer with its Intact Protein Mode and 1.8 fold increase in resolving power enables high-res analysis at high detection speed which makes it more suitable for high throughput top-down analysis.
- Aided by a new charge assignment/deconvolution algorithm, Q Exactive HF MS provides significant proteoform and protein coverage, even from a single direct infusion spectra.
- A multiplex SIM approach allows analysis of intact proteins up to about 50 kDa with isotopic resolution at LC timescale.

[www.thermoscientific.com](http://www.thermoscientific.com)

©2015 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

**Africa** +43 1 333 50 34 0  
**Australia** +61 3 9757 4300  
**Austria** +43 810 282 206  
**Belgium** +32 53 73 42 41  
**Canada** +1 800 530 8447  
**China** 800 810 5118 (free call domestic)  
400 650 5118

**Denmark** +45 70 23 62 60  
**Europe-Other** +43 1 333 50 34 0  
**Finland** +358 10 3292 200  
**France** +33 1 60 92 48 00  
**Germany** +49 6103 408 1014  
**India** +91 22 6742 9494  
**Italy** +39 02 950 591

**Japan** +81 45 453 9100  
**Korea** +82 2 3420 8600  
**Latin America** +1 561 688 8700  
**Middle East** +43 1 333 50 34 0  
**Netherlands** +31 76 579 55 55  
**New Zealand** +64 9 980 6700  
**Norway** +46 8 556 468 00

**Russia/CIS** +43 1 333 50 34 0  
**Singapore** +65 6289 1190  
**Spain** +34 914 845 965  
**Sweden** +46 8 556 468 00  
**Switzerland** +41 61 716 77 00  
**UK** +44 1442 233555  
**USA** +1 800 532 4752

**Thermo**  
SCIENTIFIC

A Thermo Fisher Scientific Brand