# Thermo Fisher

## Boosting Proteome Coverage by a Combination of the Next Generation UHPLC and a Novel Search Node in Proteome Discoverer

### Tabiwang N. Arrey<sup>1</sup>

Eugen Damoc<sup>1</sup> Bernard Delanghe<sup>1</sup>; Runsheng Zheng<sup>2</sup>; Christopher Pynn<sup>2</sup>; Oleksandr Boychenko<sup>2</sup>; Sean Woodmansey<sup>3</sup>; Wim Decrop<sup>2</sup>; Paula McBride<sup>3</sup>; Alexander Harder<sup>1</sup>

<sup>1</sup>Thermo Fisher Scientific, Bremen, Germany; <sup>2</sup>Thermo Fisher Scientific, Germering, Germany; <sup>3</sup>Thermo Fisher Scientific, Hemel Hempstead, United Kingdom

The world leader in serving science





Proteome Discoverer 3.0 software workflow



### **Abstract**

- **Thermo Fisher** SCIENTIFIC
- Due to the substantial number of MS and MS/MS scans collected during shotgun proteomics experiments, data interpretation is only possible using advanced database search engines. However, despite using advanced search algorithms, many fragment ions in the spectra remain unassigned. Some of these fragment ions may belong to peptides carrying modifications that are not considered during database search, but also to peptides that are co-isolated and co-fragmented. Here, we evaluate a novel search algorithm, CHIMERYS<sup>™</sup>, which is available in Proteome Discoverer<sup>™</sup> 3.0 software and that can identify multiple peptides in a single spectrum.

### Introduction

#### Primary Challenge

 In bottom-up proteomics MS\MS spectra often contain fragment ions from multiple co-isolated peptides, which in most cases not taken into account when doing data interpretation with existing search algorithms.

#### Prior efforts in the field

 Implementation of Iterative database search algorithms or by the use of combinations of search algorithms. However, since most of these search algorithms do not take into account that the measured intensities of fragment ions could be a sum of multiple peptides, thus, valuable information is not used during data processing. This leads to fewer peptides and protein being identification.

#### Novel approach

High-throughput methods using the Vanquish Neo UHPLC system and a PepMap Neo column in combination with a novel search algorithm, CHIMERYS to boost proteome coverage.

#### **Overview content**

- Reproducible and consistent chromatography
- Comparison of CHIMERYS to Sequest HT, Sequest HT with INFERYS rescoring for short and long gradients.

#### Summary

• CHIMERYS outperforms Sequest HT and Sequest HT algorithms with INFERYS rescoring in both semi-complex and a complex proteome for both long and short gradient. However, the highest benefits are seen in shorter gradients.

#### Trademarks

© 2021 Thermo Fisher Scientific Inc. All rights reserved. CHIMERYS™, INFERYS™ are trademarks of MSAID GmbH. SEQUEST is a trademark of the University of Washington. Powered by AWS is a trademark of Amazon, 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.

### **Materials and Methods: Sample preparation**

- Thermo Scientific<sup>™</sup> Pierce<sup>™</sup> HeLa Digest Standard (20 µg/vial) was reconstituted by adding 200 µL of 5% ACN in 0.1% formic acid (FA) in water. Sample was aspired and pipetted approximately 10 times and then transferred to an autosampler vial. 2 ul of this solution was injected on to the trapping column.
- Yeast protein Digest (Promega, 100ug) was reconstituted in 200µl of 5% ACN in 0.1% formic acid (FA) in water. It was further diluted with 0.1 % formic acid to obtain a concentration of 500 ng/ul. Sample was aspired and pipetted approximately 10 times and then transferred to an autosampler vial. 1 ul of this solution was injected on to the trapping column.

### **Materials and Methods: MS Methods**

The analytical column was connected to an Easy-Spray Source and maintained at 50 °C. Eluting peptides were sprayed on to the Thermo Scientific<sup>™</sup> Orbitrap Exploris<sup>™</sup> 480 mass spectrometer and measured.

#### 180,100 and 60 Samples Per Day, SPD

#### Trap and Elute Mode

Full Scan	
Orbitrap resolution	45,000
Scan Range	350-1,200
AGC Target (%)	300
injection time	20ms
RF Lens (%)	50
Intensity threshold	
Min. Intensity	5.00E+03
Charge state filter	
Charge state	2-5
Dynamic Exclusion	
Exclude after n times	1
Exclusion duration	25
Exclude isotopes	TRUE
Data dependent properties	
ТорМ	40
Isolation width	2
NCE (%)	28
Orbitrap Resolution	7,500
Injection time	12 ms
AGC Target (%)	50

#### 30 and 24 SPD

Irap and Elute Mode		
Full Scan		
60,000		
350-1,200		
300		
20ms		
50		
Intensity threshold		
5.00E+03		
Charge state filter		
2-5		
Dynamic Exclusion		
1		
45		
TRUE		
Data dependent properties		
30		
2		
28		
15,00		
Auto		
50		

#### 60 and 90 min gradients

#### **Direct injection**

Full Scan		
Orbitrap resolution	120,000	
Scan Range	350-1,200	
AGC Target (%)	300	
injection time	20ms	
RF Lens (%)	50	
Intensity threshold		
Min. Intensity	8.00E+03	
Charge state filter		
Charge state	2-5	
Dynamic Exclusion		
Exclude after n times	1	
Exclusion duration	45	
Exclude isotopes	TRUE	
Data dependent properties		
ТорМ	20	
Isolation width	2	
NCE (%)	28	
Orbitrap Resolution	15,000	
Injection time*	Auto	
AGC Target (%)	50	
<b>U</b> ( )		

### Materials and Methods: Data processing method

The acquired raw data files were processed with Thermo Scientific<sup>™</sup> Proteome Discoverer<sup>™</sup> 3.0 using Sequest HT, Sequest HT with INFERYS rescoring and CHIMERYS. PSM and Peptide validation is performed using Percolator.

**Thermo Fisher** 



## **Materials and Methods: LC parameters**

### LC parameters: Direct Injection mode

- Column: Thermo Scientific<sup>™</sup> Easy-Spray<sup>™</sup> PepMap<sup>™</sup> Neo 2 µm C18 75 µm X 500 mm
- Flow rate: 250 nl\min
- Gradient
  - 60 min (80 min run time)
  - 90 min (120 min run time)



60 min gradient: 8-25 % B in 50 min, 10 min to 40 %B 90 min gradient: 8-25 % B in 80 min, 10 min to 40% B

### LC parameters: Trap and Elute mode

- Easy column: Thermo Scientific<sup>™</sup> Easy-Spray<sup>™</sup> PepMap<sup>™</sup> Neo 2 μm C18 75 μm X 150 mm
- Trapping Column: Thermo Scientific<sup>™</sup> PepMap<sup>™</sup> Neo 5 μm C18 300 μm X 5 mm Trap Cartridge
- Flow rate: 1.3 to 0.4 ul/min
- Gradients:
  - 5.5 min (180 Samples Per Day, SPD),
  - 11 min (100 SPD),
  - 20.1 min (60 SPD),
  - 44.4 min (30 SPD),
  - 56.4 min (24 SPD)



Solvent A: 0.1 % FA Solvent B: 80 % ACN, 0,1%FA Weak wash solution: 0.1 % FA Strong wash solution: 80 % ACN, 0,1%FA

## **Results: Productivity and Repeatability**

 Maximizing the gradient is a key to obtaining the most out of every gradient in bottom-up proteomics. Through parallel gradient separation and sampler washing procedures along with fast sample loading on the Vanquish Neo up to 82% of the time is used to acquire useful MS spectra.



• Reproducible chromatographic and maximum MS utilization times.



## **Results: Comparison of different search algorithms**

### Yeast protein digest (500ng): Unique peptides

- Minimal variation between injections of the same gradient.
- The average number of peptides identified increased with gradient length for all the search algorithms, with CHIMERYS showing the highest identification.



### Yeast protein digest (500ng): Protein groups

• The percentage increase in the average number of protein groups identified decreases with gradient length, nevertheless CHIMERYS still outperforms the other search algorithm used in these studied.

Thermo Fishei



## **Results: Comparison of search algorithm**

### HeLa digest (200ng): Trap and Elute mode

• Also, for a complex proteome, CHIMERYS shows its superiority to the other search algorithms.



HeLa digest (200ng): Trap and Elute mode



Sequest HT INFERYS CHIMERYS

## **Results: Comparison of search algorithm**

HeLa digest (200ng): Direct Injection mode, 60 min

CHIMERYS still perform well with longer gradients



HeLa digest (200ng): Direct Injection mode, 90 min

• Even though spectra complexity is reduced by extending the gradient length, CHIMERYS is still able to confidently pick up co-fragmented precursors and identify them.



### **Conclusions**

- Due to its parallelized operation, up to 82 % MS utilization time is obtained on the Vanquish Neo with reproducible gradients and consistence results.
- Compared to Sequest HT and Sequest HT with INFERYS rescoring, the results obtained using the CHIMERYS search node show tremendous increase in the number of peptides and proteins identified from yeast- and HeLa digest.
- An increase of 30% in protein\peptide identifications for the longest gradient and of more than 80% for the shortest gradients.
- Going from 2- 4Th isolation width there observed a slight increase in the number of identified peptides.
  However, the protein group numbers did not change much.