# Improvements in LFQ for reproducible quantification of proteomic experiments: how DDA outperforms DIA

Ignacio Ortea<sup>1</sup>, Romain Huguet<sup>2</sup>, David Horn<sup>2</sup>, Andreas FR Huhmer<sup>2</sup>, and Daniel Lopez-Ferrer<sup>2</sup> <sup>1</sup> IMIBIC, Cordoba, Spain, <sup>2</sup> Thermo Fisher Scientific, San Jose, USA

# ABSTRACT

While Data Independent Acquisition (DIA) has been receiving a lot of attention lately within the proteomics community, Data Dependent Acquisition (DDA) remains the gold standard for label-free quantitation (LFQ) proteomics. DIA analyses can test whether or not a specific peptide is in a sample above a certain threshold; however, DDA methods outperform DIA when it comes to the number of peptide identifications and quantitative inter-experimental reproducibility, especially in conjugation with advanced label free quantitation software. In this work we compare HRAM quadrupole-Orbitrap <sup>™</sup> DDA, AND HRAM quadrupole-Orbitrap DIA methods head-to-head to evaluate the sensitivity and number of peptides identified and quantified ,and demonstrate that HRAM quadrupole-Orbitrap DDA technology outperforms DIA analyses significantly in proteome coverage and quantitative reproducibility.

# MATERIALS AND METHODS

#### Sample Preparation

All solvents were LC-MS grade and purchased from Fisher Scientific. Solvent A was 100% water with 0.1% formic acid. Solvent B was 80% acetonitrile, 20% water and 0.1% formic acid. Aliquots containing 500 ng/µL HELA protein digest (Pierce, PN 88328) and 1X of HRM peptide standards from Biognosys in water with 0.1% formic acid were prepared for the study.

#### LC/MS

All analyses were performed using a Thermo Scientific<sup>™</sup> EASY-nLC<sup>™</sup> 1200 system. Samples were loaded directly onto the column using the one-column (direct injection) mode, with 2µL injected onto the column, corresponding to 1 ug of total digest. The analytical columns used were a 75 µm ID Thermo Scientific<sup>™</sup> Acclaim<sup>™</sup> PepMap<sup>™</sup> column with 2 µm particles manufactured in EASY-Spray format being either 50 cm (ES803) or 75 cm in length (ES805). The column temperature was maintained at 55 °C. A linear gradient from 5% to 44 % B over 120 at 300 nL/min was used to separate the peptide mixture.

A Thermo Scientific<sup>™</sup> Q Exactive<sup>™</sup> HF MS was used. Datasets were acquired either in DDA or DIA mode.

#### Data Analysis

Raw data was processed using Thermo Scientific<sup>™</sup> Proteome Discoverer<sup>™</sup> 2.2.0.96 software. MS<sup>2</sup> spectra were searched with the SEQUEST<sup>®</sup> HT engine against a database of 42,085 human proteins including proteoforms (UniProt, May 14th, 2015). Peptides were generated from a tryptic digestion allowing for up to two missed cleavages, carbamidomethylation (+57.021 Da) of cysteine residues was set as fixed modification, and oxidation of methionine residues (+15.9949 Da), aceylation of the protein N-terminus (+42.0106) were treated as variable modifications. Precursor mass tolerance was 10 ppm and product ions were searched at 0.8 Da tolerances. Peptide spectral matches (PSM) were validated using the Percolator algorithm, based on q-values at a 1% FDR. The area of the precursor ion from the identified peptides was calculated using the new Minora Feature Detector node. Further processing was performed using the new Rt-Aligner and Feature Mapper nodes also created for the untargeted label-free quantification workflow in Proteome Discoverer 2.2. DIA data for MS1 quantitation from the Q Exactive was analyzed using Spectronaut<sup>™</sup> 9.0 software.





### RESULTS



FIGURE 3. General overview of the results window in Proteome Discoverer 2.2 software. The Peptide Group table is linked to the Consensus Features table, and the latter one is also associated to the collection of LCMS Features from each raw file. The chromatographic profiles for each LCMS Feature are shown in the Chromatogram Traces View at the bottom.

		75 cm					50 cm				
-	1	0.959	0.955	0.959	0.955	0.924	0.922	0.921	0.923	0.922	
E	0.950	1	0.96	0.961	0.961	0.932	0.932	0.929	0.932	0.93	
2 CI	0.955	0.95	1	0.95	0.957	0.927	0.926	0.923	0.927	0.925	
2	0.950	0.961	0.96	1	0.950	0.927	0.926	0.925	0.926	0.924	
	0.955	0.961	0.957	0.959	1	0.926	0.926	0.923	0.927	0.925	
	0.924	0.932	0.927	0.927	0.928	1	0.972	0.972	0.979	0.967	
ے ا	0.922	0.932	0.925	0.925	0.925	0.972	1	0.975	0.97	0.973	
0 CL	0.921	0.929	0.923	0.925	0.923	0.972	0.975	1	0.974	0.958	
2	0.923	0.932	0.827	0.926	0.927	0.979	0.97	0.974	1	0.955	
-	0.922	0.93	0.925	0.924	0.925	0.967	0.973	0.958	0.965	+	

FIGURE 5. Box-style correlation plot across the different datasets. Average correlation is ~0.96 within datasets acquired using the same column.



FIGURE 7. Scatter plot of relative protein abundance. Red dots correspond to 50 cm column and blue dots to the 75 cm one. Although it looks like the 50 cm column can provide larger dynamic range ,this result is an artifact of the roll up method, since the protein abundance is calculated after normalization as the median of the peptide abundances



**FIGURE 4.** Box plots of mean  $\log_{10}$  peptide abundance values from Protein Discoverer 2.2's LFQ algorithm for each paired comparison split out according to column length. The number of quantifiable peptides is shown on the top of each box plot.



FIGURE 6. Histogram showing the number of features used and those that have at least one missing value within the five replicates.



FIGURE 8. Venn diagram showing the number of quantified proteins that overlap between datasets, indicating good reproducibility between both separation conditions.



FIGURE 9. Histogram showing the number of precursors, peptides and protein groups identified in the DIA analyses from the library built using the Proteome Discoverer 2.2 software search results.



**FIGURE 10.** Histogram of coefficients of variation obtained from raw peptide intensities for 5 replicate datasets acquired using either DDA or DIA methods.

## CONCLUSIONS

- Data dependent acquisition in combination with a new untargeted label-free quantification workflow based on the Minora algorithm has demonstrated higher accuracy and sensitivity than data independent acquisition methods.
- The combination of the label-free quantification workflow integrated into the scaling, normalization, and study management features of Proteome Discoverer provide a powerful means for analyzing highly complex proteomics data.
- These results clearly surpass the current standards in the proteomics paradigm and rival quantitation results derived from DIA methods in terms of reproducibility and depth of analysis, but with greater efficiency, as there is no need to first generate a spectral library.
- The use of 75cm columns increased number of peptide and protein identifications, as well as the number of proteins quantified in both acquisition modes.

#### www.thermofisher.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. 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 Brazil +55 11 2730 3006 Canada +1 800 530 8447 China 800 810 5118 (ree 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 6 6885 1213 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 Sweden +46 8 556 468 00 Switzerland +41 61 716 77 00 Taiwan +886 2 8751 6655 UK/Ireland +44 1442 23355 USA +1 800 532 4752 PN10516-EN 0616S

