

Increased Dynamic Range of DDA-based label-free quantification using the CHIMERYYS algorithm

David M. Horn,¹ Kai Fritzemeier,² Katja Tham,² Frank Berg,² Daniel Hermanson,¹ Tobias Schmidt,³ Daniel Zolg,³ Martin Frejno,³ and Bernard Delanghe,² ¹Thermo Fisher Scientific, 355 River Oaks Parkway, San Jose, CA 95134, ²Thermo Fisher Scientific, Hannah-Kunath-Strasse 11, Bremen, Germany, ³MSAID, GmbH, Lichtenbergstr. 8, Garching b. München, Germany

Abstract

Purpose: Here we show how the CHIMERYYS™ intelligent search algorithm significantly increases the number of identified and quantified peptides for label-free quantification. We also show that CHIMERYYS combined with other search approaches in the Thermo Scientific™ Proteome Discoverer™ software framework can further improve results.

Methods: Human and *E. coli* proteomes were mixed and run on a Thermo Scientific™ Orbitrap Eclipse™ Tribrid™ mass spectrometer, acquiring MS/MS spectra using ion trap HCD MS/MS. Data were searched using Proteome Discoverer 3.0 software using three different processing workflows using Sequest™ HT, CHIMERYYS, and a combined 3 search node strategy, each of which also performed label-free quantification with the Minora Feature Detector.

Results: CHIMERYYS produced a significant increase in the number of identified and quantified peptides and proteins. A combined search strategy using three different search methods further improved results, producing 10,045 proteins and 144,709 unique peptides. The human and *E. coli* protein ratios were calculated with high precision.

Introduction

Current DDA-based approaches for label-free quantification match peptide identifications from a search engine to features detected independently by a quantification algorithm. With higher resolution instruments with MS1 resolution up to 480,000, it is quite common for quantification algorithms to detect substantially more than 100,000 features in a 2-hour LC/MS run. However, typical search engines can usually only identify about half of these features. Peptides may remain unidentified due to unexpected post-translational modifications of enzyme cleavage specificity, but quite frequently peptides are simply not identified because they are co-isolated and fragmented with several other peptides to produce a chimeric spectrum.

The CHIMERYYS intelligent search algorithm helps address this issue by identifying significantly more peptides in chimeric spectra, dramatically increasing the number of peptide identifications in complex proteomics LC-MS/MS runs. For high resolution MS/MS data, CHIMERYYS can identify more than 10 PSMs in the same MS/MS spectrum. We show here that the pairing of CHIMERYYS with the Minora feature detection algorithm in Proteome Discoverer 3.0 software leads to an increase in the number of quantified peptides and proteins in complex proteomics data compared to a Sequest HT-based approach. We further show that CHIMERYYS can be combined with other search approaches including INFERYYS™ Rescoring to further increase the number of identified and quantified peptides.

Materials and methods

Increasing amounts of an *E. coli* tryptic digest sample (0, 25, 50, 100, and 200 ng) were mixed with a human tryptic digest to make a total of 1 µg sample per mixture. Each sample was run in triplicate using an Orbitrap Eclipse Tribrid mass spectrometer equipped with a Thermo Scientific™ FAIMS Pro™ interface, using a 2-hour gradient and two compensation voltages (-55 and -70 V). MS scans were detected at 240K resolution and HCD MS/MS scans were collected in the ion trap using a 0.7-*m/z* isolation window with a 1 s cycle time.

All data analysis was performed using Proteome Discoverer 3.0 software. The first analysis used a Sequest HT-based Processing workflow for label-free quantification (Figure 1). The Spectrum Files RC node performs mass recalibration for both the precursors masses for the Sequest HT search and the LCMS Features created by the Minora Feature Detector. The nodes used a combination of the March 2022 canonical human and *E. coli* proteome FASTA databases downloaded using the Proteome Discoverer annotation service.

Data were also analyzed using a CHIMERYYS-based Processing workflow (Figure 2). The CHIMERYYS workflow includes a "Top N Peaks Filter" node prior to the search, passing through the top 20 peaks per 100 *m/z* window for MS/MS spectra. The filter significantly improves CHIMERYYS results for ion trap MS/MS data but is not recommended for high resolution MS/MS data. The CHIMERYYS search node used a fragment tolerance of 0.3 Da, fixed Carbamidomethyl cysteine and oxidized methionine, and the inferyys_2_1_fragmentation prediction model.

A standard label-free quantification Consensus workflow was used for both workflows using default parameters except requiring a channel occupancy of 50% in the Precursor Ions Quantifier node. *p*-values were calculated using the background-based method and ratios were calculated using the pairwise ratio approach. All protein, peptide group, and PSM counts are reported at a 1% FDR threshold, using Percolator for PSM-level FDR for all 3 search approaches, qvality for peptide group FDR, and the Protein FDR Validator node for Protein FDR.

Figure 1. Sequest HT-based processing workflow for label-free quantification.

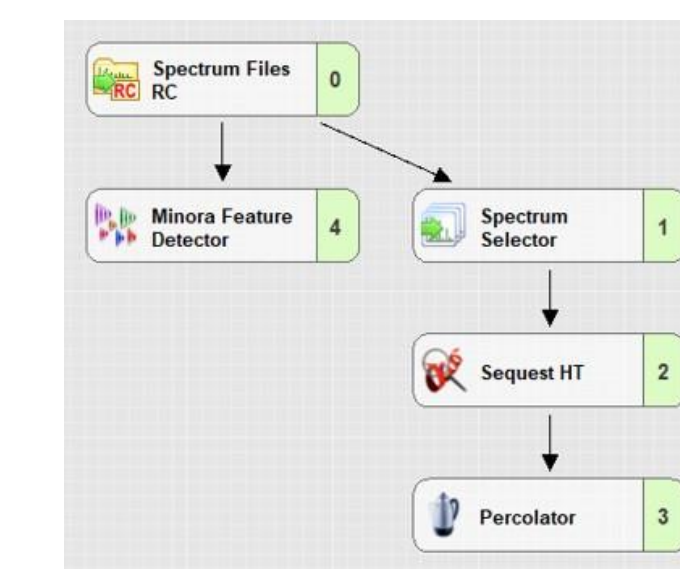
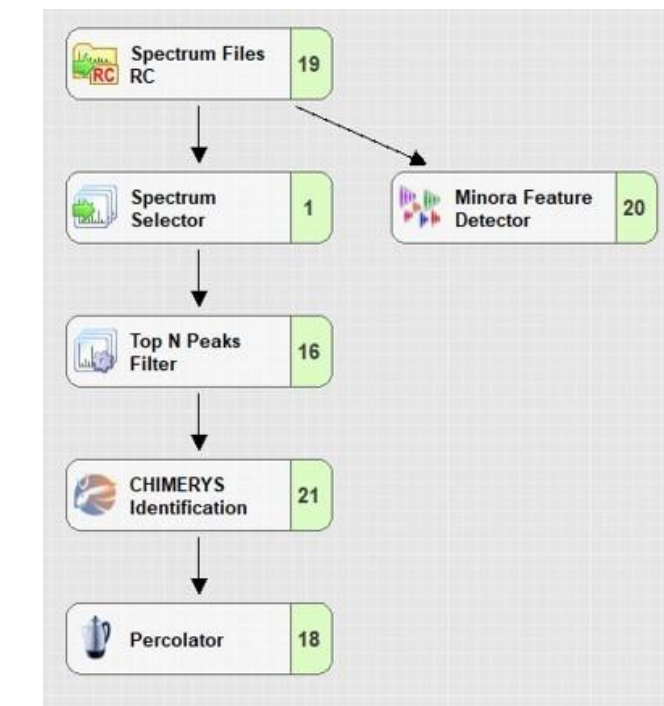
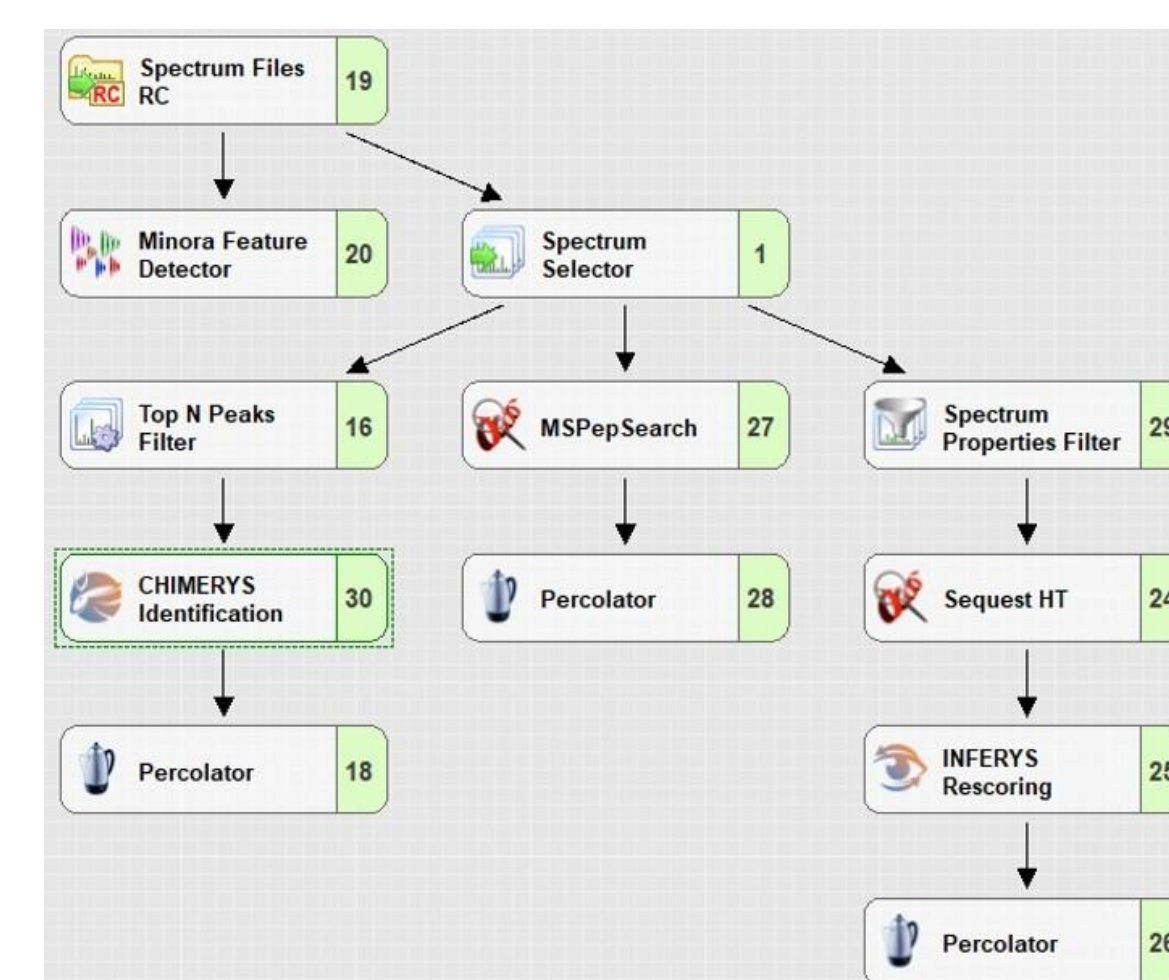


Figure 2. CHIMERYYS-based processing workflow for label-free quantification.



Data were subsequently analyzed using a branched workflow. While CHIMERYYS currently supports enzyme-specific cleavage and peptides with fixed carbamidomethyl cysteine, TMT™, or TMTpro modifications, other nodes can be used to identify peptides produced by semi-specific cleavage and peptides with post-translational modifications. Sequest HT with INFERYYS™ Rescoring¹ has been shown to significantly increase the number of peptide identifications for peptides produced by non-specific cleavage. To speed up the search, the FASTA databases for the human and *E. coli* proteomes were pre-indexed prior to search. Another branch of the workflow included the MSPepSearch algorithm using the NIST OT HCD library to identify proteins commonly detected in human samples. The combined processing workflow with all 3 search nodes is shown in Figure 3.

Figure 3. Branched processing workflow for label-free quantification with 3 search nodes and the Minora Feature Detector. The Sequest HT and INFERYYS Rescoring search were used to identify semi-tryptic peptides for both the human and *E. coli* proteomes, while the MSPepSearch node was used to identify commonly detected peptides with post-translational modifications.



Results

Figure 4. Comparison of proteins, unique peptides, and PSMs between the Sequest HT and CHIMERYYS searches.

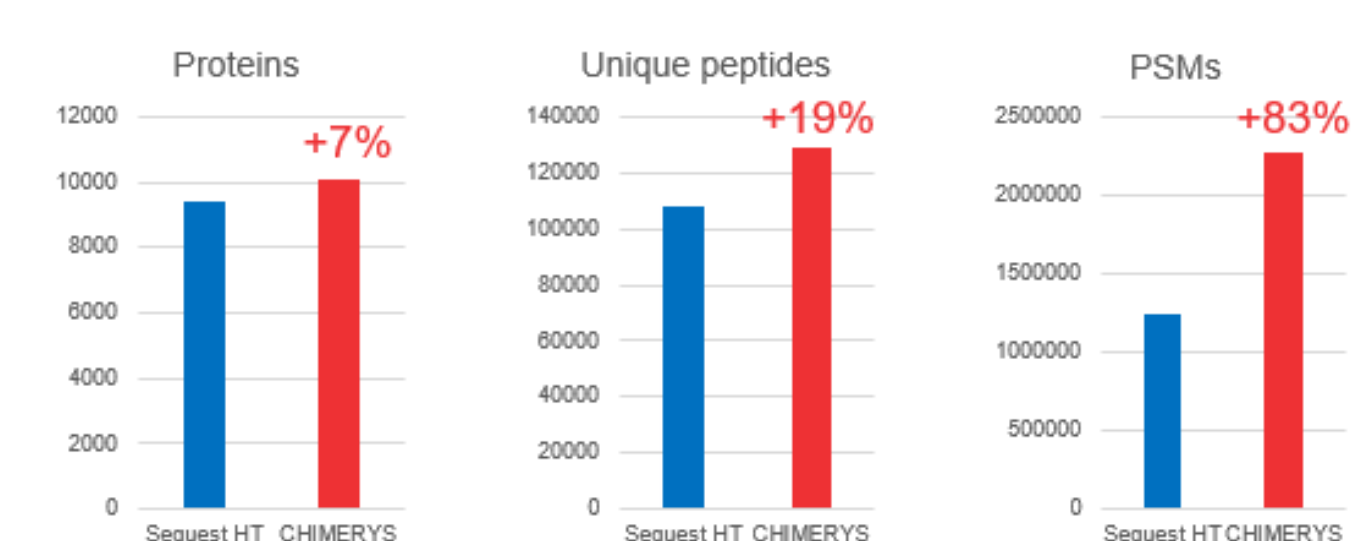
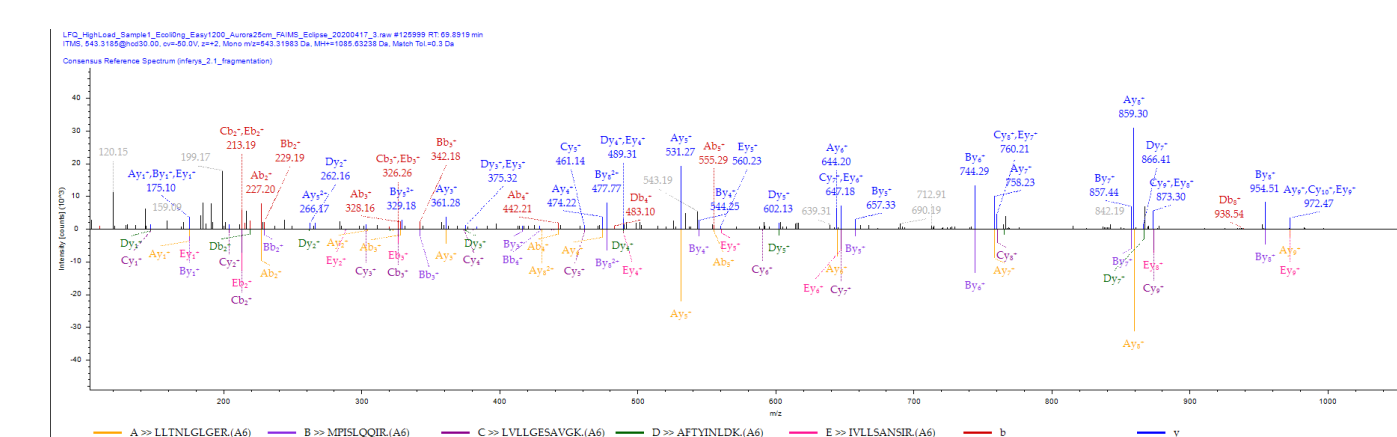


Figure 4 shows a comparison of the protein and peptide identification results from the Sequest HT and CHIMERYYS workflows. CHIMERYYS produces 19% more unique peptides and 7% more proteins compared to Sequest HT, with an equal increase in protein identifications for both the human and *E. coli* proteomes. While this increase is less than the 50-70% observed for high resolution MS/MS data, the improvement is still significant given the nominal resolution and mass accuracy of ion trap MS/MS data. The CHIMERYYS search produced up to 6 PSMs in a single MS/MS spectrum while the Sequest HT search can only identify 1 PSM. Figure 5 shows one example of an MS/MS spectrum where 5 peptides were simultaneously identified by CHIMERYYS.

Figure 5. Fragment Match Spectrum for an MS/MS scan with 5 PSMs identified using CHIMERYYS. The lower spectrum in the mirror plot shows the INFERYYS-predicted combined MS/MS spectrum for all 5 peptides. Almost every abundant peak in the spectrum has been assigned as a fragment.



The number of quantified *E. coli* proteins by search engine is shown in Figure 6. For each concentration level, the CHIMERYYS results produce about the same 7% increase in the number of quantified proteins as is seen for the number of protein IDs.

Figure 6. Number of quantified *E. coli* proteins by amount of sample. There are about 6% more quantified proteins using CHIMERYYS at each concentration level. The number of quantified human proteins are similar across all samples and, thus, not displayed here. CHIMERYYS quantified about 7% more human proteins for each sample compared to Sequest HT.

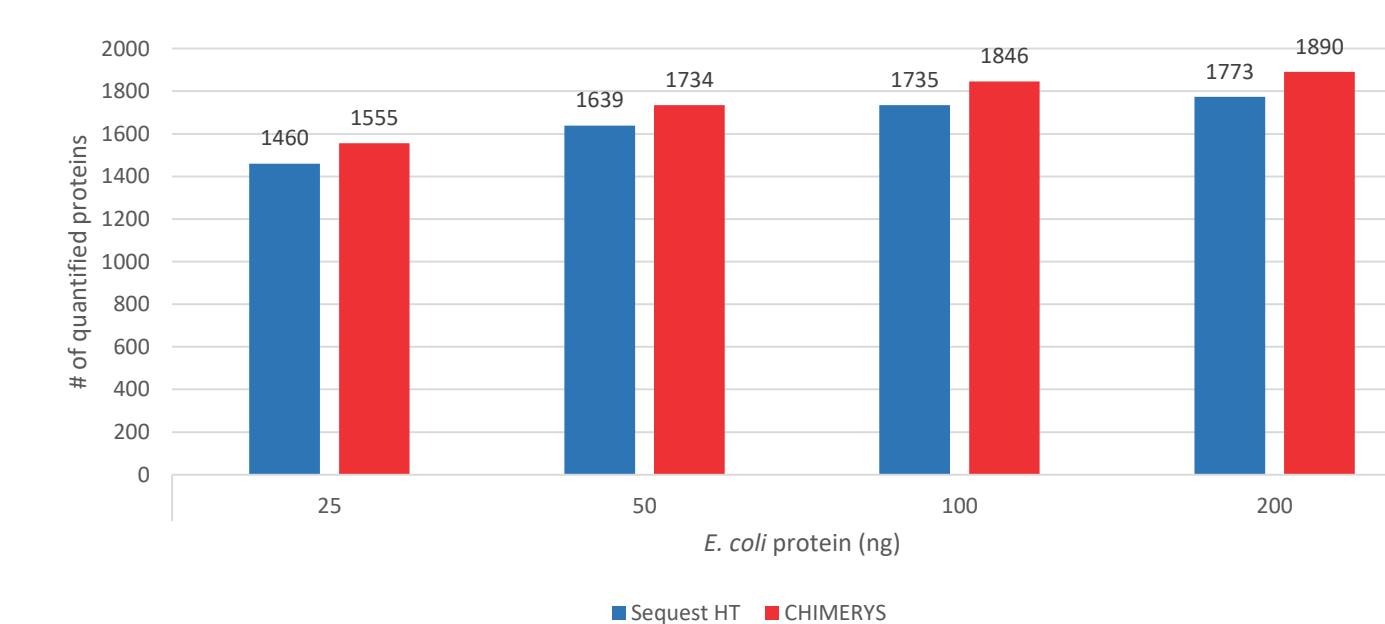
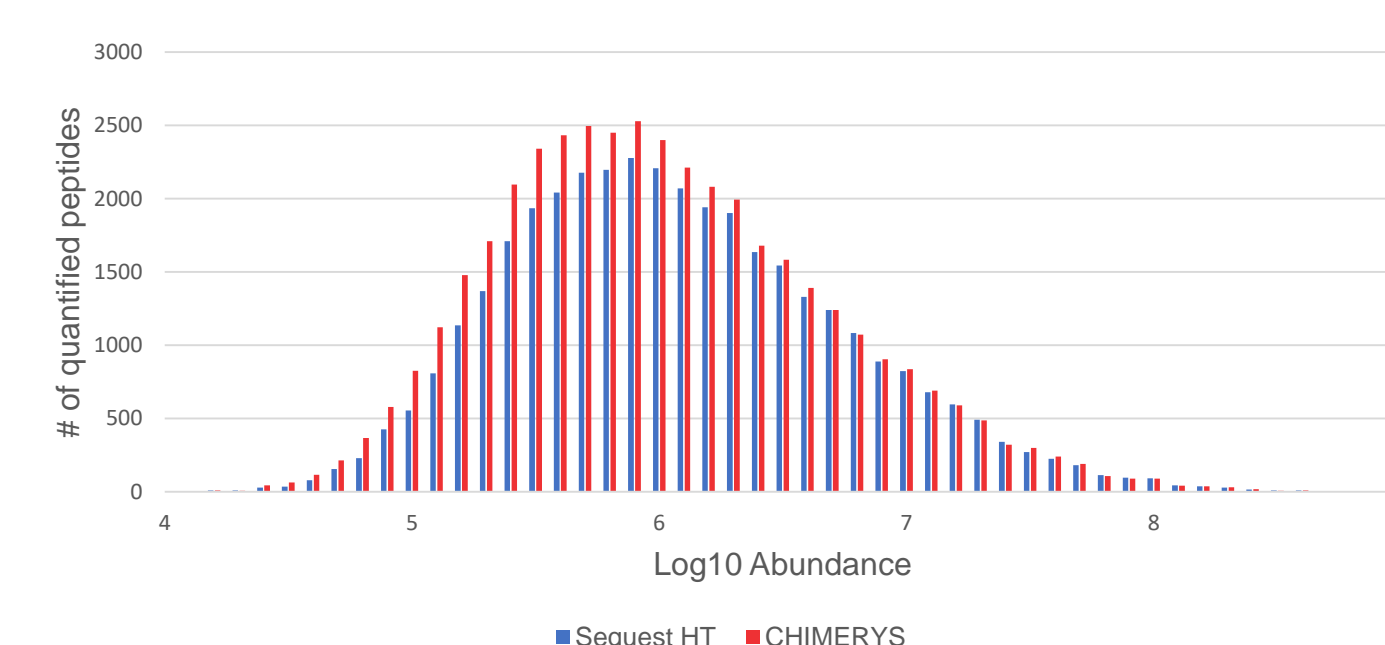


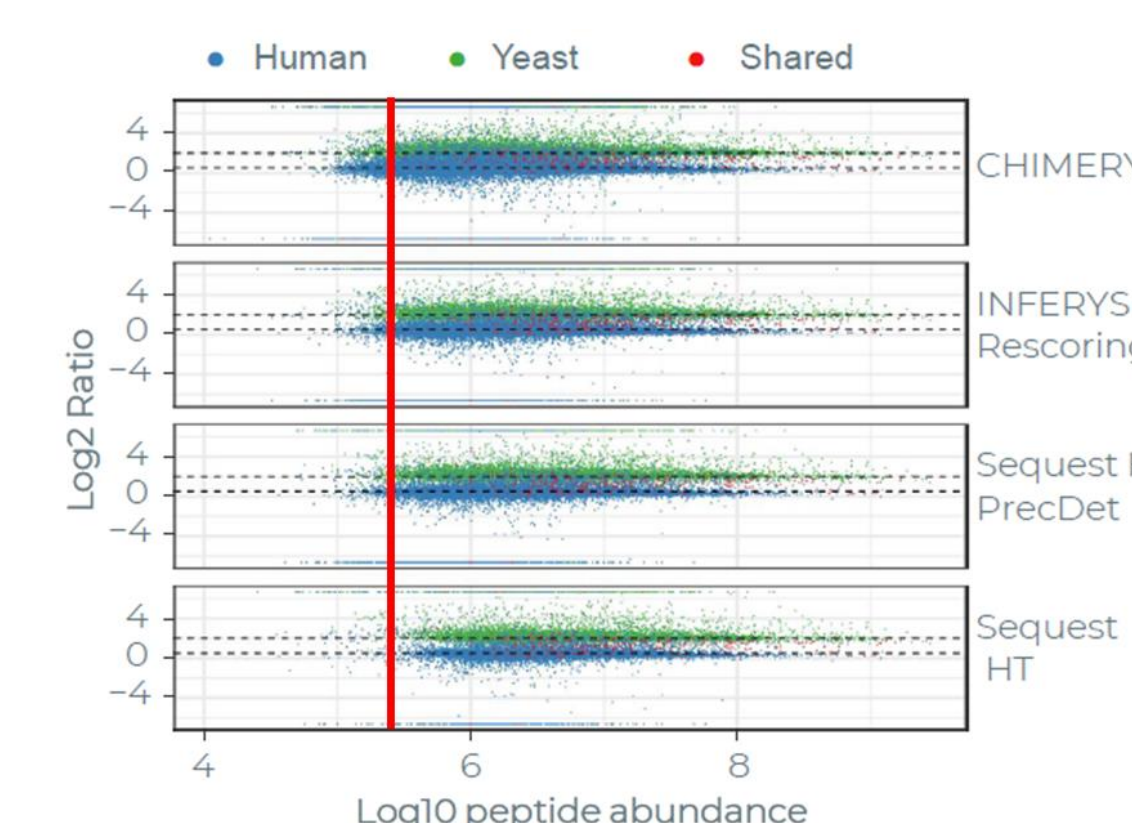
Figure 7 shows the distribution of *E. coli* peptide abundances for the three 200 ng *E. coli* datasets. The higher abundance peptides are quantified by both workflows, while the lower abundance peptides are more frequently quantified using the CHIMERYYS-based workflow.

Figure 7. Distribution of peptide group abundances for the 200 ng *E. coli* samples. CHIMERYYS identifies more low abundance peptides with Log10 intensities below 6.



The dynamic range is improved due to the ability of CHIMERYYS to identify lower abundance peptides in mixed spectra. For nominal mass, this improvement is relatively subtle, but can be seen more clearly when searching high resolution MS/MS data. Figure 8 shows the improvement in the dynamic range for identification using a high-resolution MS/MS dataset with human and yeast proteomes. The CHIMERYYS results show the high dynamic range of quantification.

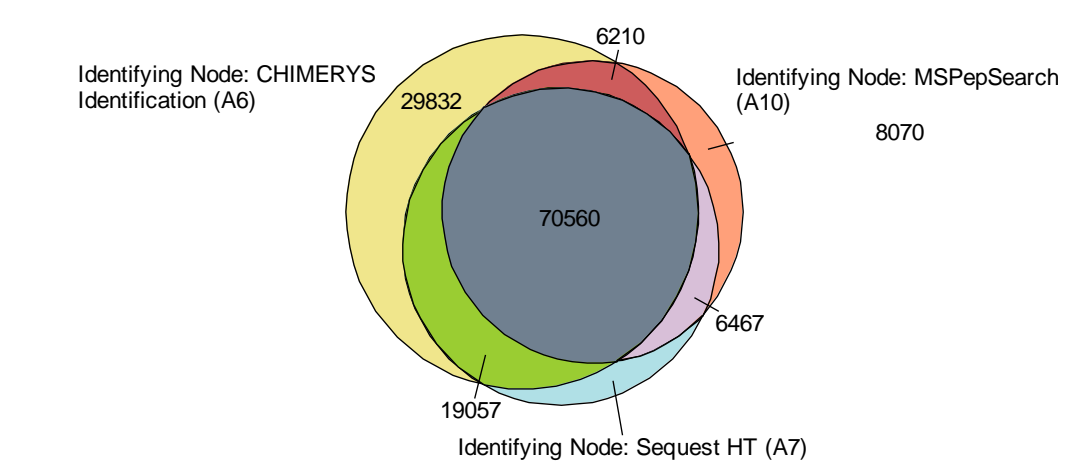
Figure 8. Distribution of peptide group abundances for a human/yeast proteome mixture acquired using Orbitrap-based HCD MS/MS and processed with different search strategies.



Results from combined 3 search node method

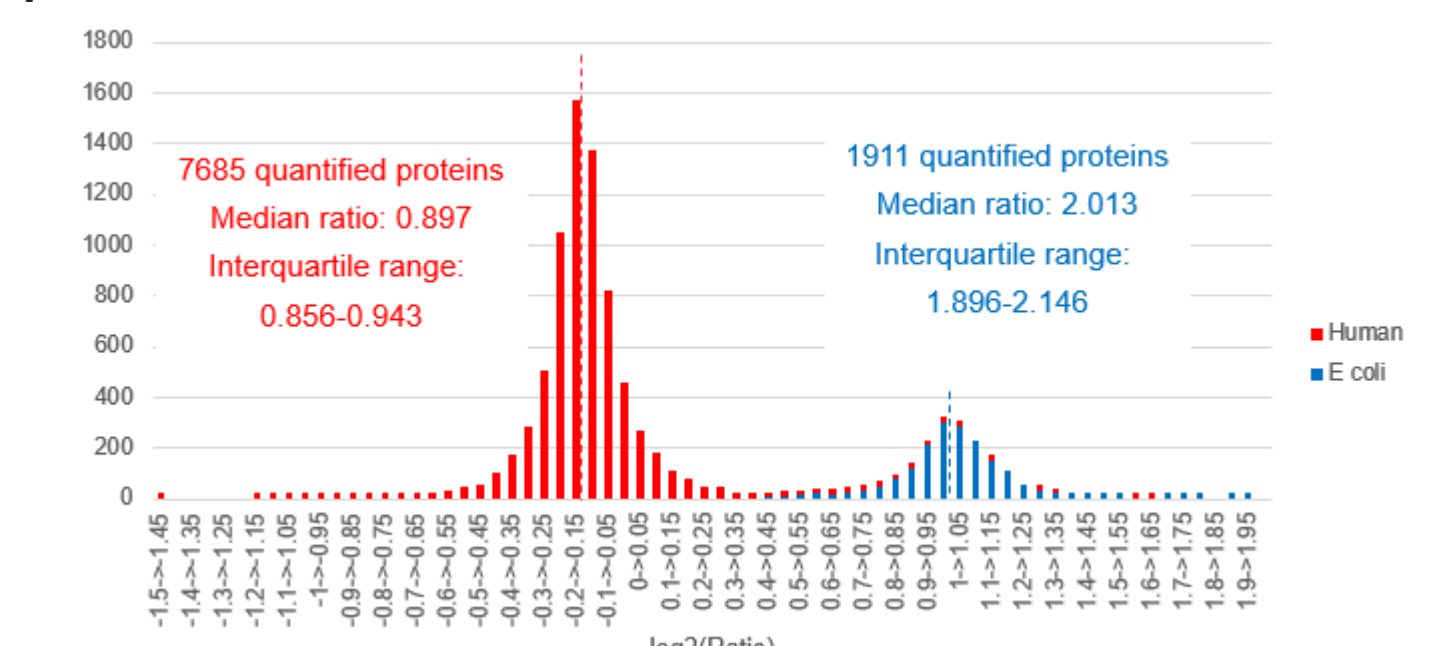
The combined workflow produced a similar number of protein identifications (10045) as the CHIMERYYS search alone (10120), but there was a moderate increase in the number of unique peptide identifications (+12%). The overlap in identifications between the three different search strategies is shown in Figure 9. Many of the identifications that are unique to Sequest HT and INFERYYS Rescoring correspond to semi-tryptic peptides and many unique to MSPepSearch results include modified peptides.

Figure 9. Venn diagram showing the overlap in unique peptide IDs for the 3 different search nodes. While CHIMERYYS produced by far the most identifications, both MSPepSearch and Sequest HT identified several thousand peptides not identified by CHIMERYYS.



The Minora Feature Detector and Feature Mapper nodes were improved for the Proteome Discoverer 3.0 software release to better support CHIMERYYS search results. As a result, there is substantially reduced false quantification in Proteome Discoverer 3.0 software than past releases for label-free quantification. Figure 10 shows the distribution of quantification values for the human and *E. coli* proteins for the comparison of the 200 ng and 100 ng *E. coli* samples.

Figure 10. Histogram of log2 ratios for the human and *E. coli* proteins for the 3 search node results. The dashed lines show expected ratios.



Of the 144700 identified peptides, 104916 had measured abundances for the 200 ng samples, 106738 had measured abundances for the 100 ng samples, 106193 had measured abundances for the 50 ng samples, 104112 for the 25 ng samples, and 97247 for the 0 ng samples. The lower abundance samples are expected to have fewer quantified peptides due to the decreasing amounts of *E. coli* peptides.

Conclusions

CHIMERYYS produces a substantial increase in identifications compared to a standard Sequest HT search. CHIMERYYS is not only able to identify more peptides in these mixed spectra, but also the additional peptides in the same spectra are usually lower in abundance than the peptide identified by Sequest HT. Thus, when CHIMERYYS is paired with the Minora Feature Detector, the workflow will be able to identify lower abundance peptides than a traditional search engine. Subsequent work after these data were acquired shows that CHIMERYYS can identify even more peptides from ion trap MS/MS data of human samples using a wider isolation window (1.5-*m/z*) and a higher maxIT (15 ms) than what was used for these data (0.7-*m/z* and 10 ms). We would thus expect another substantial increase in identified and quantified peptides using this updated wide window acquisition method that takes better advantage of the CHIMERYYS algorithm.

References

- "INFERYYS rescoring: Boosting peptide identifications and scoring confidence or database search results", Rapid Commun. Mass Spectrom., 2021

Trademarks/licensing

© 2022 Thermo Fisher Scientific Inc. All rights reserved. CHIMERYYS and INFERYYS are trademarks of MSAID, GmbH. SEQUEST is a trademark of the University of Washington. TMT is a trademark of Proteome Sciences plc. 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.