New peak detection (Pyco) and isotope grouping (Prism) algorithms for an improved compound detection workflow

too low.

Yovany Cordero Hernandez, Pedro Navarro, Konstantin Ayzikov, Martin Strohalm, James Dillon, Alexander Tiegel, Ute Comberg, Ralf Tautenhahn, Alexander Makarov, Christoph Henrich, Thermo Fisher Scientific, 355 River Oaks Parkway, San Jose, CA, 95134

Abstract

A reliable peak detection algorithm is key for any compound detection workflow. Here we introduce the new detection workflow in Thermo Scientific[™] Compound Discoverer[™] 3.3 software, based on a novel parameter-free peak detection algorithm, Pyco, and a redesigned isotope grouping algorithm, Prism. We show that this new approach works more consistently on various chromatography methods, as peak integration does not rely on restrictive peak model assumptions, such as Gaussian shape or symmetry. Using peak quality metrics, we corroborate that both identification and quantification results have been significantly improved. We were able to reduce false positive compounds by avoiding over-splitting of peaks. The new detection workflow is not only more precise, but also more efficient, reducing the overall computational time for processing.

Trace quality filtering

Gaps in traces are interpolated when gaps are found between two non-zero values and the gap is no longer than two missing points.





Data Reduction Process



Introduction

Challenges in peak detection

- Overlapping peaks/deconvolution
- Baseline detection and removal
- Accurate peak integration
- Peak-model-based detection and -integration vs. model-free

Baseline separated peaks



Problems with model-based peak integration



- Sensitivity vs. Specificity
- "Real" peak or just noise ?
- How to measure peak quality?

Baseline removal using AirPLS



Trace Denoising and Smoothing



Isotope Detection & Grouping

Prism (Pattern Recognition of Isotopes in Small Molecules)

- Fast isotope pattern detection tuned for small molecule data.
- Looks for isotopes by mass shifts relative to an A_0 mass.
- Supports detection of the following isotopic elements: carbon, hydrogen, oxygen, nitrogen, sulfur, bromine, chlorine



Results

Improved sensitivity and runtime

- New peak detection is faster
- New peak detection is more sensitive: detects more compounds and more ions per compound
- Result tables can be much larger \rightarrow filter based on Peak Rating

| Compound Discoverer Software version | Intensity threshold | Peak Rating Threshold | Number of Compounds after Group Compounds, <u>unfiltered</u> | Number of Compounds after Group Compounds, <u>filtered by Peak</u> <u>Rating</u> in Compound Discoverer Software | Runtime Detect Compounds |
|---|------------------------|--|--|---|----------------------------------|
| 3.2 | 10,000 | n/a | 59,958 | Not possible to filter by Peak Rating | 17h 3min |
| 3.3 | 10,000 | Peak Rating ≥ 4 in at least 3 samples | 104,332 | 19,931 | 4h 27min |
| | | | | | |

121 samples (soy sauce). Processed on a HP Z840, 2 x (Intel Xeon CPU E5-2667 @3.2 GHz, 8 cores), 64GB RAM, purchased in 2017.

High baseline, noisy peak

Noisy peak, low signal

Partially overlapping peaks



Compound Detection workflow

Compound Detection in Compound Discoverer software is performed in four steps.

- Trace Generation and -filtering a.
- Detection of chromatographic peaks b.
- Detection of isotopic patterns C.
- Grouping of adducts and fragments d.





Peak Detection

Peak Detection algorithm in images







Isotope detection using PRISM



Step 2. Load mass spectrum at chromatographic peak apex. 266.0113 367.1505 421.0612 00 200 300 400 500 600

Select window around the A_0 mass. Step 4. Look for <u>carbon isotopes</u> that match

the mass offset and expected intensity of current A₀ mass.

Step 5. Look for non-carbon isotopes that Non-C Isotopes match the mass offset and expected intensity of the current A_0 mass.

Step 3.

150.9692 z=0

157.9743 z=1

164.9485 z=0



Compare list of detected isotopes to list of detected chromatographic peaks. Select matching chromatographic peaks and group with the A_0 peak.

Quantification Reproducibility

- Evaluation using data published by Li et al.³, 1100 spiked compounds at known concentrations
- Publication contains comparison of free software tools and a much older version of **Compound Discoverer** software (v2.1, current version is 3.3) CVs calculated with

compounds detected in

at least 2 raw files



Legacy (CD 3.2)

Pyco (CD3.3)

- Pyco shows a significantly reduced variability compared to the legacy algorithm in Compound Discoverer software
- Pyco's variability is on par with the best results from other software compared in the paper

Conclusions

- Increased sensitivity reduces chances of missing compounds
- Peak Rating is a flexible mechanism to filter out low quality compounds and adds an additional metric for evaluation
- New peak detection contributes to improved quantification
- New detection workflow improves workflow run time

References

1.Z.-M. Zhang, S. Chen, and Y.-Z. Liang, Baseline correction using adaptive iteratively reweighted penalized least squares.

Measures the first unexpected change in direction of intensity to detect splitting and integration of multiple peaks.

Final Peak Rating

Captures shape quality by

measuring the normalized

its neighbors on either side.

variance between a point and

Captures shape quality by

calculating the number of

of intensity vector.

Peak Rating values in Compound Discoverer software are calculated based on the four different Peak Quality Factors, but also take into account CV and relative peak area.

Adduct Grouping

Analyst 135 (5), 1138-1146 (2010).

2. Adapted from *Pandey et al.*, Metabolomics (2020), MetaClean: a machine learning-based classifier for reduced false positive peak detection in untargeted LC–MS metabolomics data

3. Li et al., Anal Chim Acta. (2018), Comprehensive evaluation of untargeted metabolomics data processing software in feature detection, quantification and discriminating marker selection.

Trademarks/licensing

© 2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.