

A High-Throughput Integrated HRAM-MS Method Enables IGF-1 Quantification, Targeted Variants Monitoring, and Untargeted Variants Screening in a Single Injection

Yu Zhou, Xiaolei Xie, Kristine, Van Natta, Bradley Hart, Shen Luan, Debadeep Bhattacharyya
Thermo Fisher Scientific, 355 River Oaks Parkway, San Jose, CA 95134

ABSTRACT

Purpose: To develop a fast, robust, and highly integrated HRAM data acquisition and data analysis workflow for serum IGF-1 on a modified Thermo Scientific™ Orbitrap Exploris™ mass spectrometer. With fast scanning and enhanced mass accuracy using EasyIC™, this novel method enables IGF-1 quantification, targeted variants identification, and unknown variants screening in a single injection. In addition, the integrated data analysis provides complementary information for improving data quality and confidence.

Methods: We utilized a two-step protein precipitation, combined with online sample extraction on a Thermo Scientific™ Vanquish™ Duo UHPLC. Analytes were detected on a modified orbitrap mass spectrometer. In a single method, data is acquired using a) multiplexed SIM scan mode for IGF-1 quantification, b) tMS² mode for targeted known variants confirmation, and c) full scan mode for unknown variants screening.

Results: The modified orbitrap shows good analytical sensitivity, linear dynamic range and excellent reproducibility for reliable serum IGF-1 measurements. Intact IGF-1 was quantified at a LLOQ of 10 ng/mL with a standard curve range of 15.6–2,000 ng/mL ($r^2 > 0.99$).

INTRODUCTION

Insulin-like growth factor 1 (IGF-1) quantification has many clinical research and anti-doping applications. Differentiating IGF-1 variants is important for capturing mutations and correct interpretation of false results. High-Resolution Accurate-Mass (HRAM) MS methods have unique advantages for reliably measuring and differentiating IGF-1 species, attributes which might be missed by immunoassays and non-HRAM methods. Using the most abundant m/z values in the $[M+7H]^+$ isotopic cluster as a quantifier ion, QTOF based HRAM methods have been successfully developed for IGF-1 quantification. However, this has practical challenges in a true routine context for differentiating IGF-1 and its variants due to limited mass resolving capability, compromised specificity and quantitative performance (using ± 10 ppm extraction windows), and mass deviation issues.

MATERIALS AND METHODS

Sample Preparation

Calibrators were prepared in 20% acetonitrile in 1X phosphate buffered saline with 1% BSA. Internal standard N₁₅-labeled IGF-1 was purchased from Prospect International.

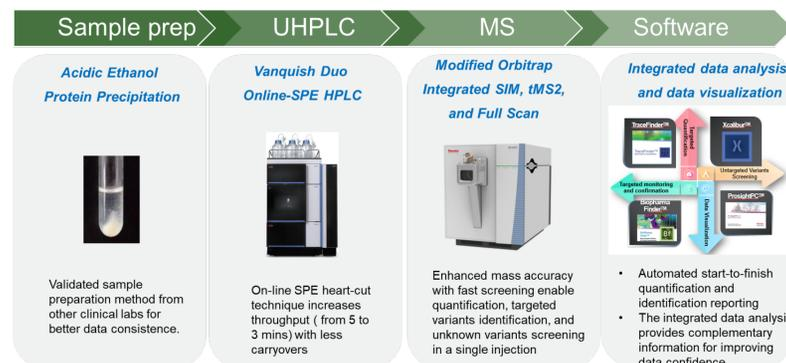
Test Method(s)

Internal standard (IGF-1-N¹⁵) were added to the serum samples which were first precipitated using acidic ethanol, and second cryoprecipitated after neutralization. Automated online SPE cleanup and chromatographic separation was performed on a Thermo Scientific™ Vanquish™ Duo UHPLC, followed by analyte detection on a modified Thermo Scientific™ Orbitrap Exploris™ mass spectrometer. A Thermo Scientific™ Hypersil Gold™ C8, 50 x 2.1 mm, 5 μ m column was used for separation. Total run time was 3 minutes.

Data Analysis

Automated quantification of SIM data was performed by Thermo Scientific™ TraceFinder™ software. Top down and intact protein analysis were processed using Thermo Scientific™ BioPharma Finder™ and ProSight™ software.

Figure 1. Enhanced mass accuracy and fast scanning enable high-throughput integrated IGF-1 quantitation and its variants identification in serum



RESULTS

Figure 2. Mass extraction (± 2.5 ppm extraction window) of a single isotope from IGF-1 in the 7+ charge state (m/z 1093.5225) was used to generate extracted ion chromatograms for accurate quantification

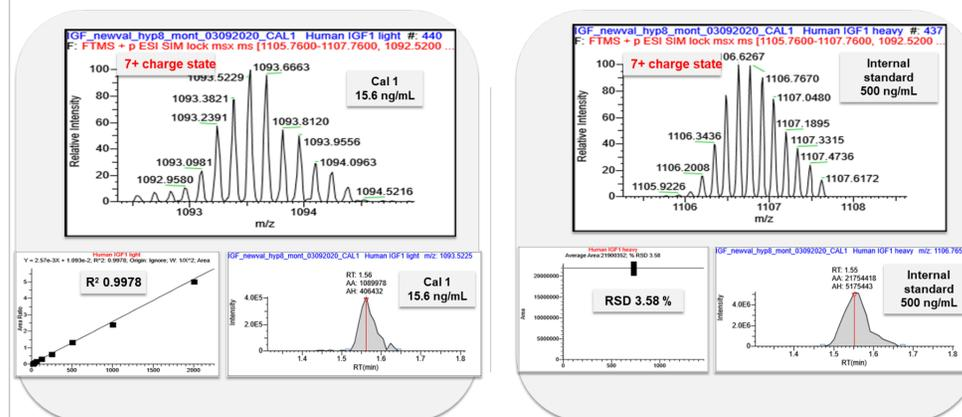


Figure 3. EASY-IC enable IGF-1 mass accuracy within ± 0.8 ppm in Calibrators and ± 1 ppm in serum samples across 130 hours of non-stop run time.

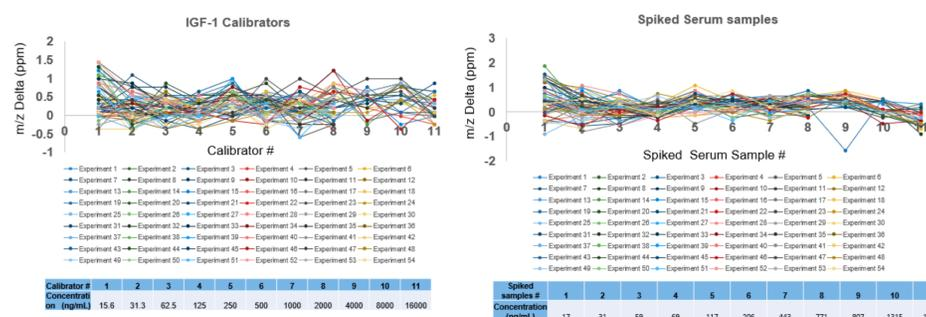


Figure 4. Overlay of 54 Calibration Experiments across 130 hours of non-stop run time for the IGF-1 quantification robustness.

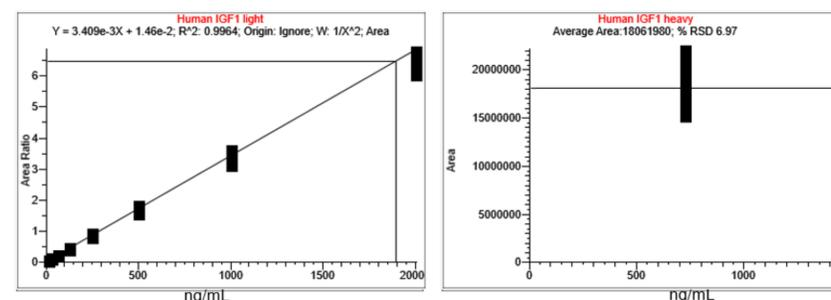


Table 1. Robustness statistics for replicates (n=54) of a calibration curve across 130 hours of non-stop run time for IGF-1.

Level (n=54)	Concentration	% Con Bias (Average)	% Ion Ratio	% CV
Cal1	15.6 ng/mL	-2%	pass	7%
Cal2	31.3 ng/mL	2%	pass	4%
Cal3	62.5 ng/mL	2%	pass	4%
Cal4	125 ng/mL	2%	pass	3%
Cal5	250 ng/mL	2%	pass	4%
Cal6	500 ng/mL	1%	pass	3%
Cal7	1000 ng/mL	-1%	pass	3%
Cal8	2000 ng/mL	-6%	pass	3%

Table 2. Robustness statistics for replicates (n=54) of 7 spiked serum samples across 130 hours non-stop run time for IGF-1.

Spiked sample # (n=54)	Concentration	% Recovery (Average)	% Ion Ratio	% CV
Sample 1	15.6 ng/mL	-4%	pass	13%
Sample 2	31.3 ng/mL	-5%	pass	6%
Sample 3	62.5 ng/mL	-7%	pass	4%
Sample 4	125 ng/mL	-7%	pass	3%
Sample 5	250 ng/mL	-16%	pass	3%
Sample 6	500 ng/mL	-12%	pass	3%
Sample 7	1000 ng/mL	-20%	pass	3%

CONCLUSIONS

- The Vanquish Duo UHPLC platform combined with the modified Orbitrap mass spectrometer provides a fast, robust, and highly integrated HRAM data acquisition and data analysis workflow for serum IGF-1.
- Normal-type, known variants and unknown IGF-1 variants can be measured in a single run.
- Based on accurate mass and signature isotope distribution, HRAM full scan mode allows identification of IGF-1 and its variants, including pathogenic, mutations, and post translational modifications. A great advantage to a full scan monitoring is the ability to re-analyze data when unknown mutations are suspected.
- The Vanquish Duo provided both samples clean-up and analytical separation in a 3-minute method, enhancing the ability of the mass spectrometer to detect compounds.
- The modified orbitrap mass spectrometer provides excellent robustness of data over long run-times without maintenance.

REFERENCES

- Maus, A., Kemp, J., Milosevic, D., Renuse, S., Pandey, A., Singh, R. J., Grebe, S. K. G., Center of Mass Calculation in Combination with MS/MS Allows Robust Identification of Single Amino Acid Polymorphisms in Clinical Measurements of Insulin-Like Growth Factor-1. *J. Proteome Res.* 2020, 19, 186-193 DOI: 10.1021/acs.jproteome.9b00494.
- Grübner M., Greco G., Flexible HPLC instrument setups for double usage as one heart-cut-2D-LC system or two independent 1D-LC systems, Thermo Fisher Scientific Technical Note 73298

TRADEMARKS/LICENSING

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All 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.