

Assessing Biosimilarity by Monitoring Multiple Critical Quality Attributes of an Intact Monoclonal Antibody Drug Using Orbitrap Native LC-MS

Aaron O. Bailey¹, Bobby Gamez¹, Terry Zhang¹, Stephane Houel¹, Shanhua Lin², Guanghui Han³, Katie Southwick¹, Wendy Sandoval³, Jonathan L. Josephs¹. ¹Thermo Fisher Scientific, 355 River Oaks Parkway, San Jose, CA, USA, 95134; ²Thermo Fisher Scientific, 1228 Titan Way, Sunnyvale, CA, USA, 94085; ³Department of Microchemistry, Proteomics and Lipidomics, Genentech Inc., South San Francisco, CA, USA, 94080

ABSTRACT

Purpose: To demonstrate feasibility of workflow for monitoring multiple critical quality attributes of a therapeutic protein drug as analyzed in the intact form.

Methods: We analyzed proteins using pH-elution ion exchange coupled directly to native ESI-MS. Isoform identification was accomplished using Thermo Scientific™ BioPharma Finder™ software and XIC-based isoform relative abundance measurements were performed using Thermo Scientific™ Chromeleon™ software.

Results: We demonstrate that ion exchange native MS provides high precision intact mass measurement of Herceptin® and a trastuzumab biosimilar and that Chromeleon software may be used to measure the XIC-based relative abundances of individual CQA isoforms.

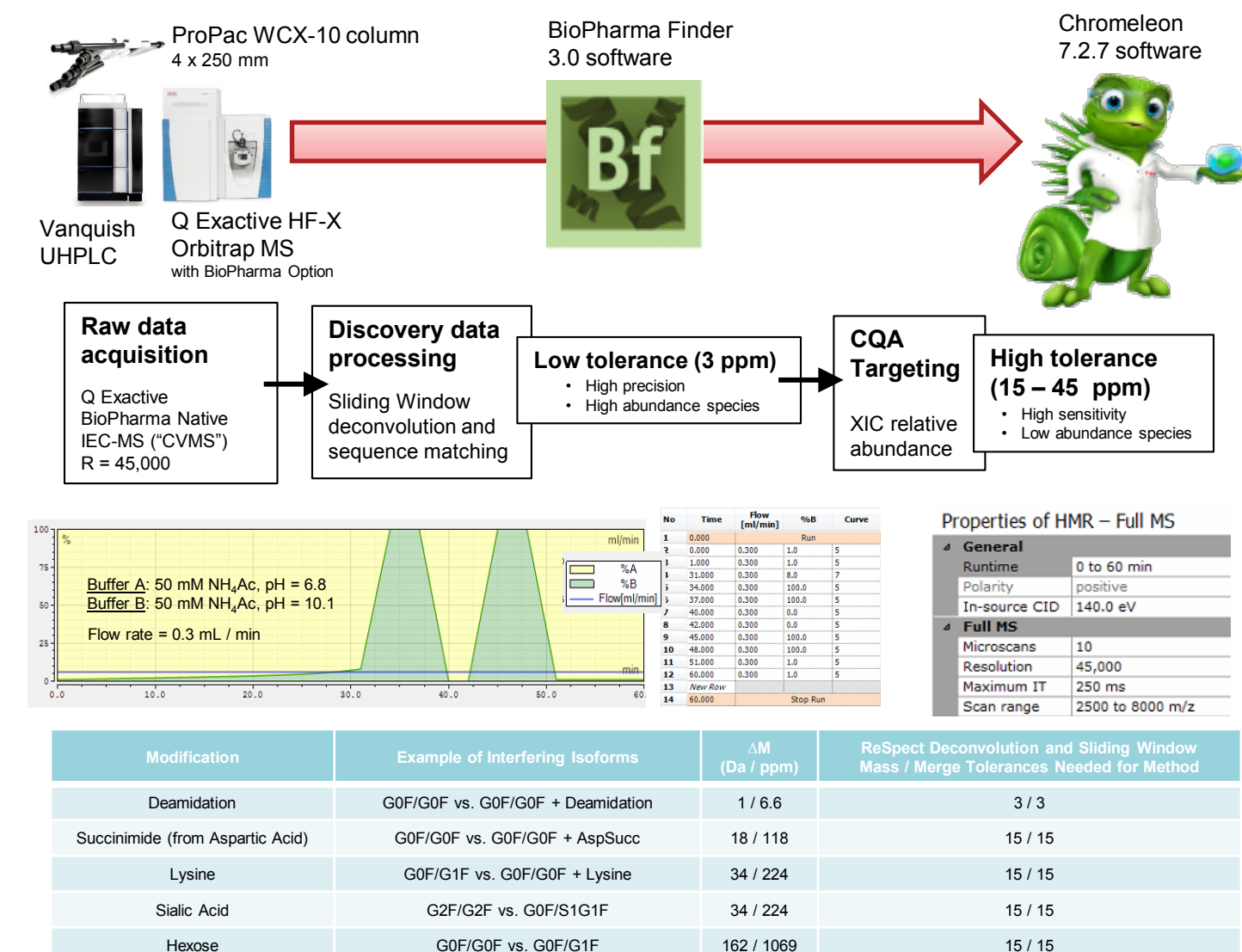
INTRODUCTION

Manufacturing of innovator biologics can be successfully mimicked to produce generic “biosimilar” drug products. In order to satisfy safety and efficacy requirements, a biosimilar drug must be reasonably comparable to an innovator. Comparability is directly assessed by measuring a panel of critical quality attributes (CQAs). Recently, a multi-attribute method has been demonstrated to measure several CQAs simultaneously using LC-MS peptide mapping data. Additional CQAs, such as charge variants or size variants (i.e., aggregation/fragmentation), are conventionally monitored at the intact level using ion exchange (IEC) or size exclusion chromatography (SEC), respectively. Native LC-MS approaches, such as SEC-MS and IEC-MS, combine the traditional gold standard chromatographic assays with accurate mass measurement to allow isoform-specific monitoring of multiple CQAs of intact therapeutic proteins.

MATERIALS AND METHODS

Intact trastuzumab was separated by IEC (Thermo Scientific™ ProPac™ WCX-10 column) directly coupled to MS. Mobile phases consisted of aqueous 50 mM ammonium acetate, using a pH gradient from 6.8 to 10.1. LC-MS was accomplished using a Thermo Scientific™ Vanquish™ H-Class UHPLC system with a variable wavelength detector directly connected to a Thermo Scientific™ Q Exactive™ HF-X Orbitrap™ mass spectrometer. Native LC-MS raw data were analyzed using a time-resolved deconvolution approach utilizing Sliding Window and ReSpect algorithms in BioPharma Finder software. Chromeleon software was used to integrate XIC peak areas and relatively compare CQAs of innovator and biosimilar drug samples.

Figure 1. Platform for on-line native IEC-MS method. Essential instrumentation for our platform includes the ProPac WCX-10 weak cation exchange column, the Vanquish Horizon UHPLC, the Q Exactive HF-X Orbitrap mass spectrometer, and BioPharma Finder data analysis software.

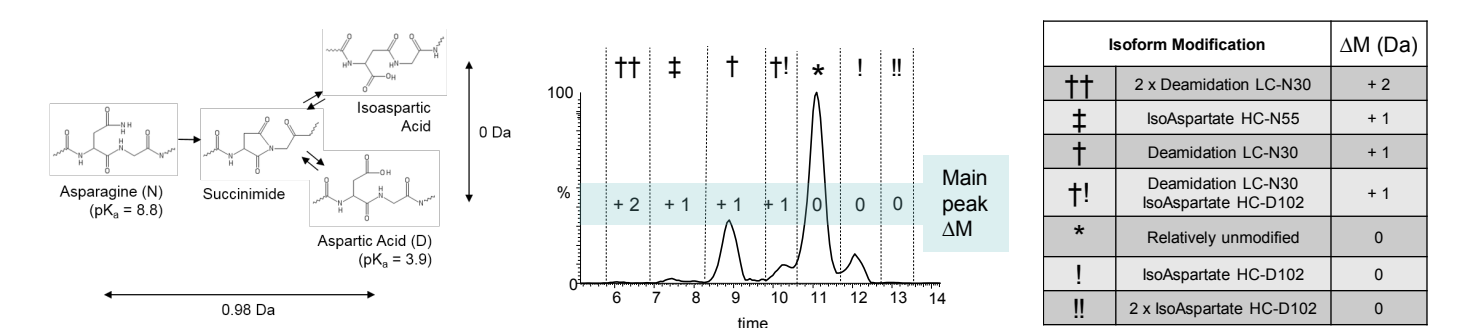


DEAMIDATION / ISOMERIZATION PATHWAY DOMINATES TRASTUZUMAB CHARGE VARIANT MS PROFILE

Deamidation and isomerization isoforms pose challenge of near-isobar heterogeneity

Trastuzumab includes several known residues in the light chain and heavy chain sequences, which may undergo deamidation (LC-N30) or isomerization (HC-D102) or both at once (HC-N55). These modified isoforms are considered to be “near-isobars”, which exhibit mass differences in the range of 0-2 Da, and pose an issue for accurate mass identification by MS. These deamidation and isomerization isoforms are observed in the charge variant profile and can be observed directly using pH-elution ion exchange coupled directly to Orbitrap native MS.

Figure 2. Asparagine deamidation and aspartic acid isomerization are observed as distinctly separated chromatographic peaks in the charge variant profile. These isoforms are near-isobars, ranging in theoretical mass differences of 0-2 Da, making independent intact mass analysis a challenging endeavor. Isoform identities were previously attained by Harris et al., using fractionation and peptide mapping of each LC peak (Reference 1).



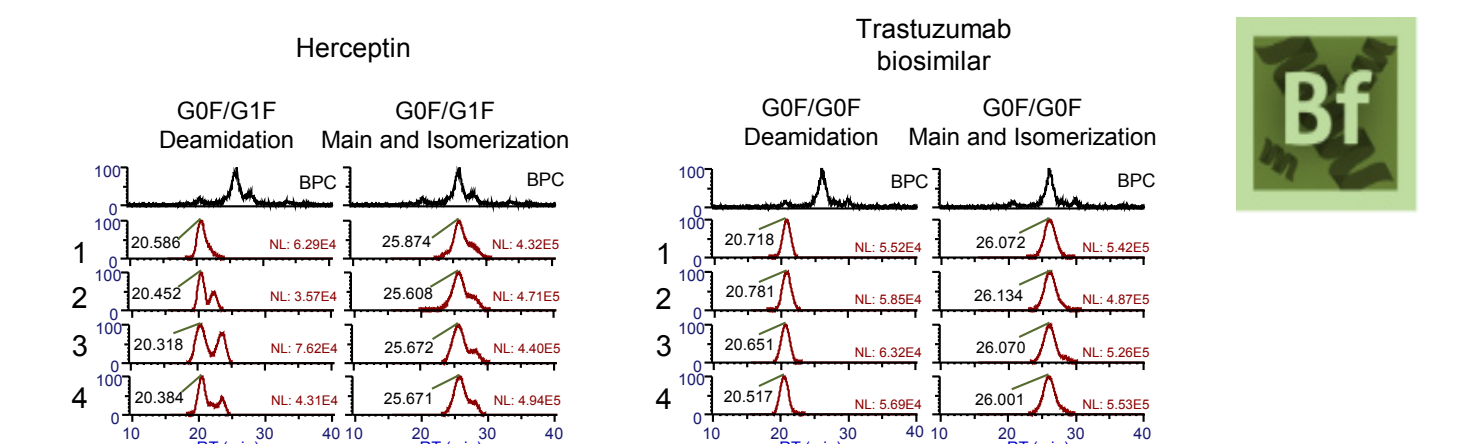
LOW TOLERANCE SLIDING WINDOW DECONVOLUTION (3 ppm) ALLOWS SUB-Da PRECISION FOR INTACT MASS MEASUREMENT

High precision of Orbitrap mass analyzer is demonstrated by reproducible results.

We injected 1 μ L (20 μ g) volumes of trastuzumab samples in formulation buffer and acquired native LC-MS data on a Q Exactive HF-X system. The results were analyzed using ReSpect and Sliding Window algorithms in BioPharma Finder software. We found a reproducible difference (4 replicates per sample) of ~1 Da between the abundant glycoforms corresponding to the acidic (deamidation) and main (relatively unmodified) LC peaks. Interestingly, we also observed a further ~1 Da difference between innovator Herceptin and a trastuzumab biosimilar, which we could not immediately explain.

Figure 3. High precision measurements show that an approximately 1 Da mass shift is reproducibly measured for abundant glycoforms at the main vs. acidic peaks, which is highly consistent with deamidation.

Sample	Modification (isoform)	Mean Average Mass (Da)	Avg. Mass Standard Dev. (Da) (ppm)	Mean Mass Accuracy vs. Theoretical (Da) (ppm)	Mean Relative Abundance (%)	Mean Apex RT (min)
Herceptin	2xAG2GF	148057.95	0.28 / 1.85	3.06 / 20.23	63.93	25.94
Herceptin	1xDeamidation (N), 2xAG2GF	148058.36	0.83 / 5.46	3.50 / 23.09	4.76	21.34
Herceptin	1xAG2GF, 1xAG2GF1F	148219.81	0.18 / 1.19	2.86 / 18.89	100.00	25.71
Herceptin	1xDeamidation (N), 1xAG2GF, 1xAG2GF1F	148221.33	0.43 / 2.85	3.41 / 22.48	10.63	20.44
Herceptin	2xAG2GF	148381.81	0.39 / 2.59	2.79 / 18.44	70.36	25.49
Herceptin	1xDeamidation (N), 2xAG2GF1F	148383.05	0.99 / 6.55	3.05 / 20.15	5.08	20.58
Trastuzumab Biosimilar 1	2xAG2GF	148058.74	0.15 / 1.01	3.87 / 25.55	100.00	26.07
Trastuzumab Biosimilar 1	1xDeamidation (N), 2xAG2GF	148060.26	0.30 / 1.95	4.42 / 29.16	7.87	20.67
Trastuzumab Biosimilar 1	1xAG2GF, 1xAG2GF1F	148220.65	0.19 / 1.25	3.71 / 24.50	40.16	25.89
Trastuzumab Biosimilar 1	1xDeamidation (N), 1xAG2GF, 1xAG2GF1F	148222.84	0.65 / 4.26	4.95 / 32.67	4.44	20.57
Trastuzumab Biosimilar 1	2xAG2GF1F	148382.58	0.37 / 2.44	3.58 / 23.62	15.76	25.74
Trastuzumab Biosimilar 1	1xDeamidation (N), 2xAG2GF1F	148383.60	0.91 / 6.00	3.62 / 23.87	0.90	20.32



MODERATE TOLERANCE SLIDING WINDOW DECONVOLUTION (15 ppm) ALLOWS SENSITIVE AND CONFIDENT DETECTION OF LOW LEVEL CQA ISOFORMS

Confident detection of low-level charge variant CQAs on intact mAb

Native IEC-MS analysis of intact trastuzumab resulted in a chromatographic profile consistent with previously published data. Time-resolved deconvolution analysis of the innovator drug resulted in identification of several specific isoforms comprised of differential N-glycosylation in combination with a variety of low level charge-imparting PTMs, such as deamidation, C-terminal lysine removal, or sialylation of N-glycans. We confirmed that the main peak consisted of several isoforms corresponding to variable N-glycosylations of the trastuzumab amino acid sequence. When comparing the innovator and biosimilar products, we found marked differences in glycoform profile measured at the main peak. Furthermore we found measurable differences in several charge variant isoforms when comparing innovator and biosimilar. These differences observed were consistent with a comparison of these same two samples via MAM HRAM peptide mapping analysis (data not shown).

Figure 4. Glycoform profiles are easily identified using ReSpect and Sliding Window algorithms in BioPharma Finder software. Our “moderate” mass tolerance of 15 ppm is wide enough to detect both deamidated and main peak forms of trastuzumab. We readily observed distinct glycoform profiles for Herceptin and biosimilar trastuzumab samples.

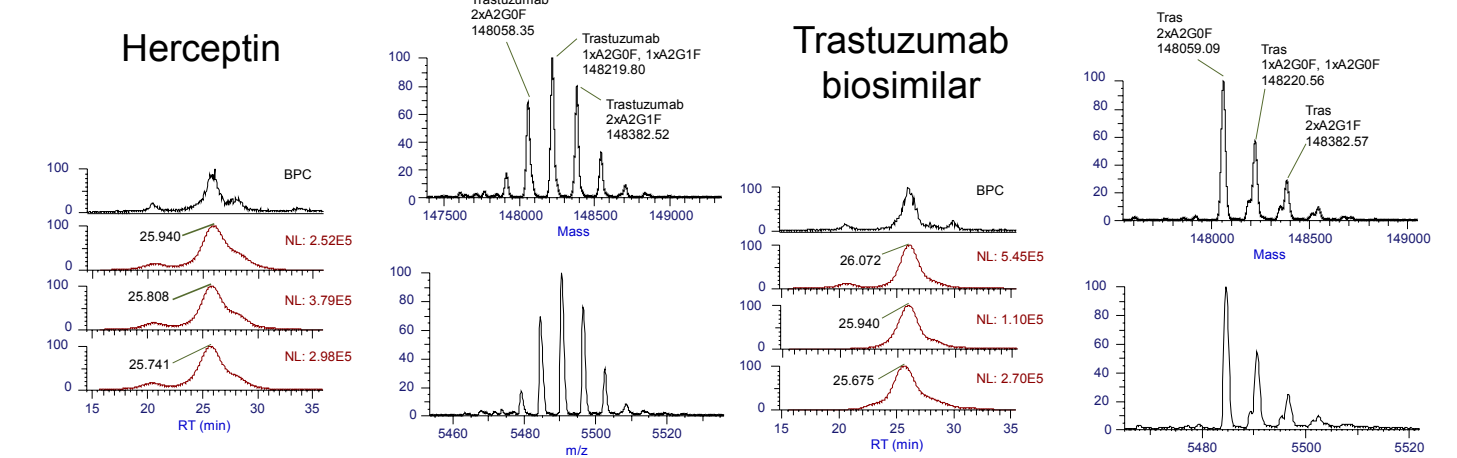
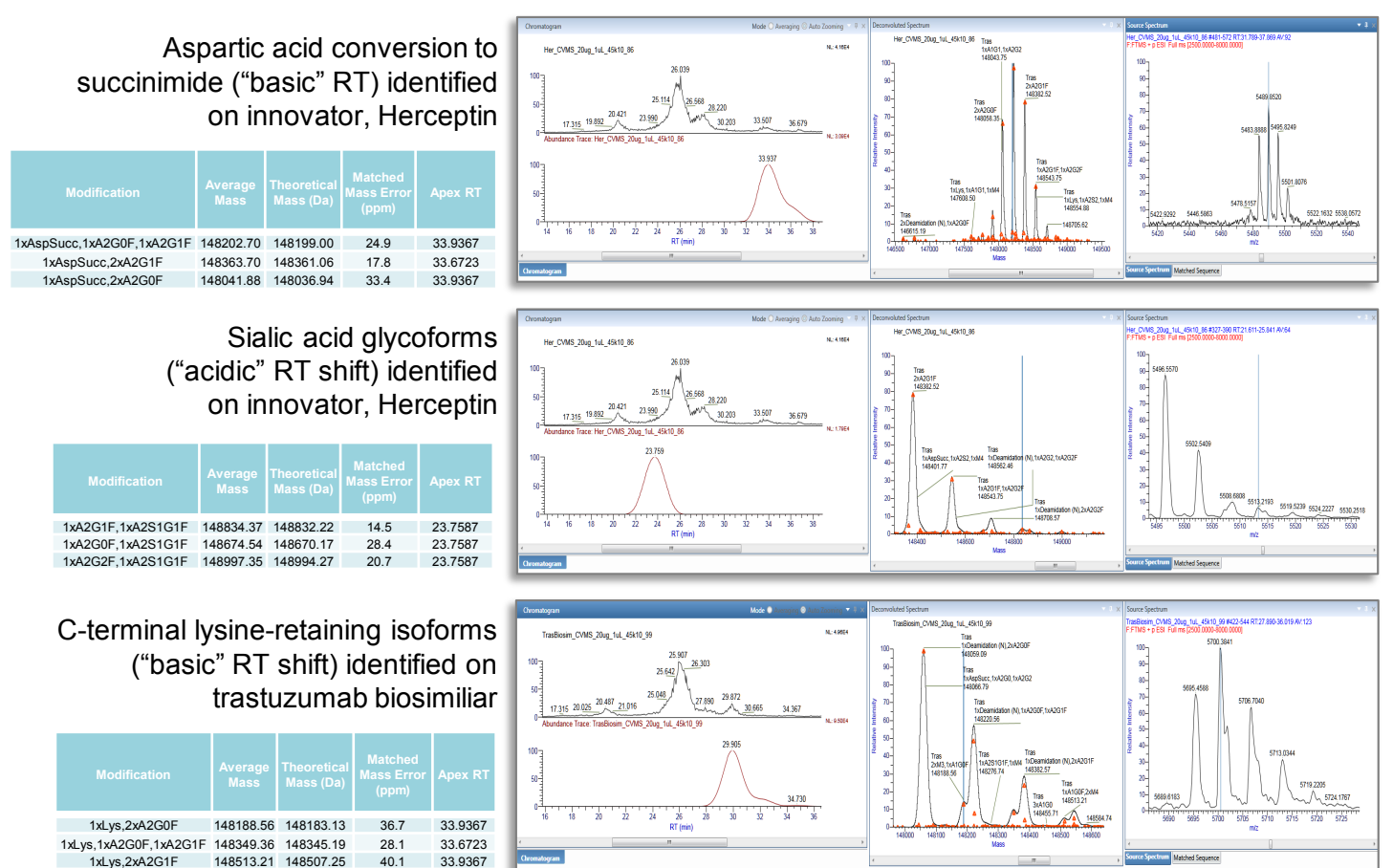


Figure 5. Screenshots from BioPharma Finder software show that lower-level CQA isoforms such as succinimide conversion from aspartic acid, sialic acid glycoforms, and C-terminal lysine-retaining isoforms are distinctly identified by mass and retention time shift relative to the main peak.



MONITORING CQA ISOFORMS IN CHROMELEON SOFTWARE

Large scale relative abundance measurements of isoforms identified by BioPharma Finder software

Quantitative monitoring of protein drugs can be easily accomplished using peptide mapping LC-MS using a so-called Multi-Attribute Method (MAM) (Reference 2). This type of approach allows CQAs to be become characterized and then monitored in a routine fashion. The combination of BioPharma Finder software and Chromeleon software serves as a powerful framework for MAM applications to enable deep protein characterization and further method streamlining for downstream process development.

We sought to develop a workflow that performs targeted quantitation of isoform charge states identified by Sliding Window deconvolution of native IEC-MS datasets. We manually exported the centroid m/z values of the charge states of isoforms identified by ReSpect / Sliding Window analysis in BioPharma Finder software. Centroid m/z values were averaged from the 4 replicates and these average values were targeted in Chromeleon.

Figure 6. BioPharma Finder intact protein analysis results can be curated to determine which species should be monitored as CQAs. Interesting isoforms can be selected (box checked) to mark for export in spreadsheet format. To export the constituent charge states’ centroid m/z values for the isoforms identified by BioPharma Finder software, right-click > “Export Checked” > “All Levels”.

Level	Charge State	Intensity	m/z Centroid	Set as Reference Component
1	Charge State	25	3.82E+05	Export All
2	Charge State	28	6.84E+05	Export Checked
3	Charge State	27	7.13E+05	Component Level Only
4	Charge State	28	3.48E+05	All Levels
5	Charge State	29	1.71E+05	All Levels

Figure 7. Isoform charge state m/z values and RT information can be used to create a processing method with flexible options for XIC integration. We quantified 3-4 charge states for each isoform, which is shown here in an overlay plot shaded for customized integration specific for the main peak and deamidated isoforms of G0F/G0F, extracted at 3, 15, 30, and 45 ppm tolerances.

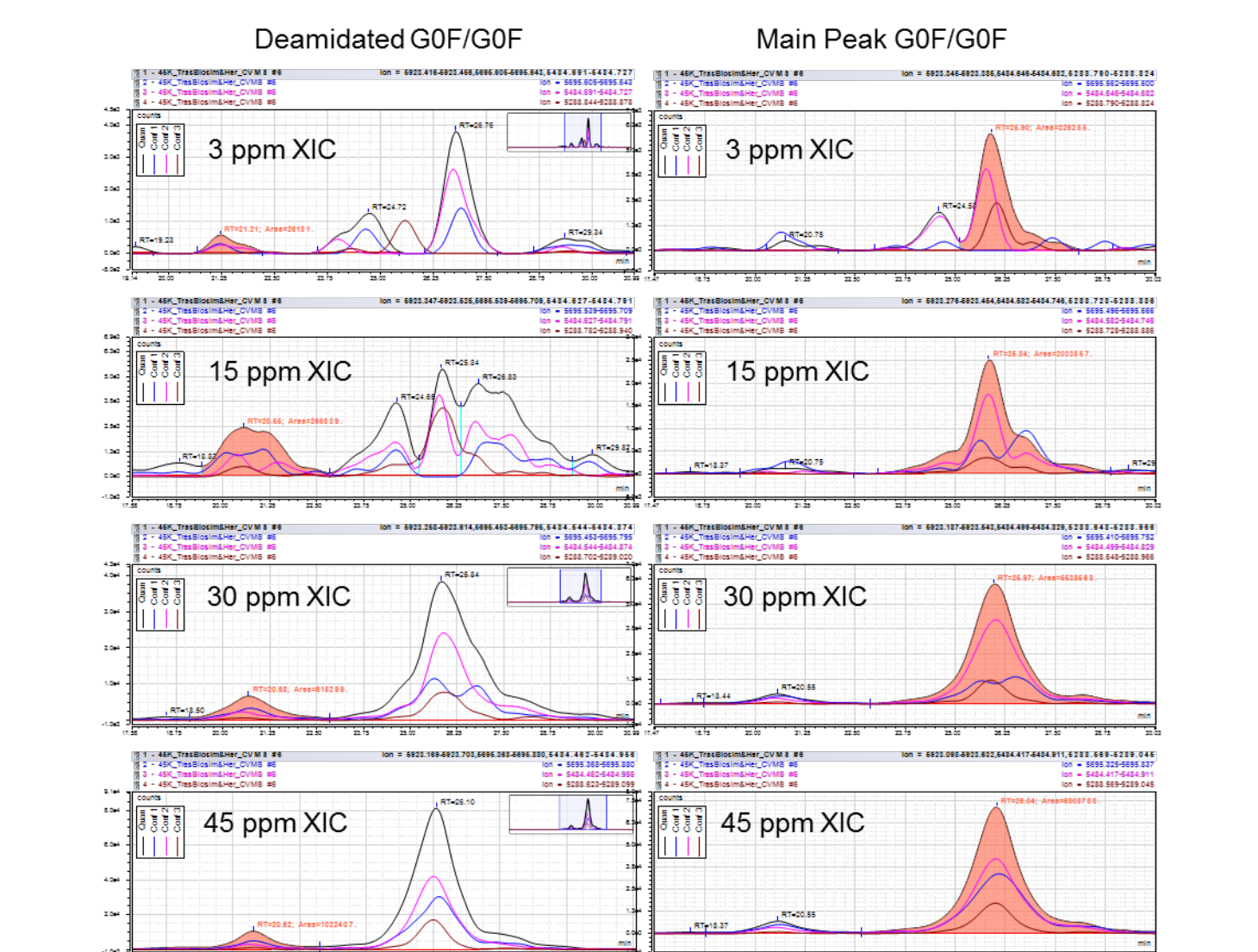
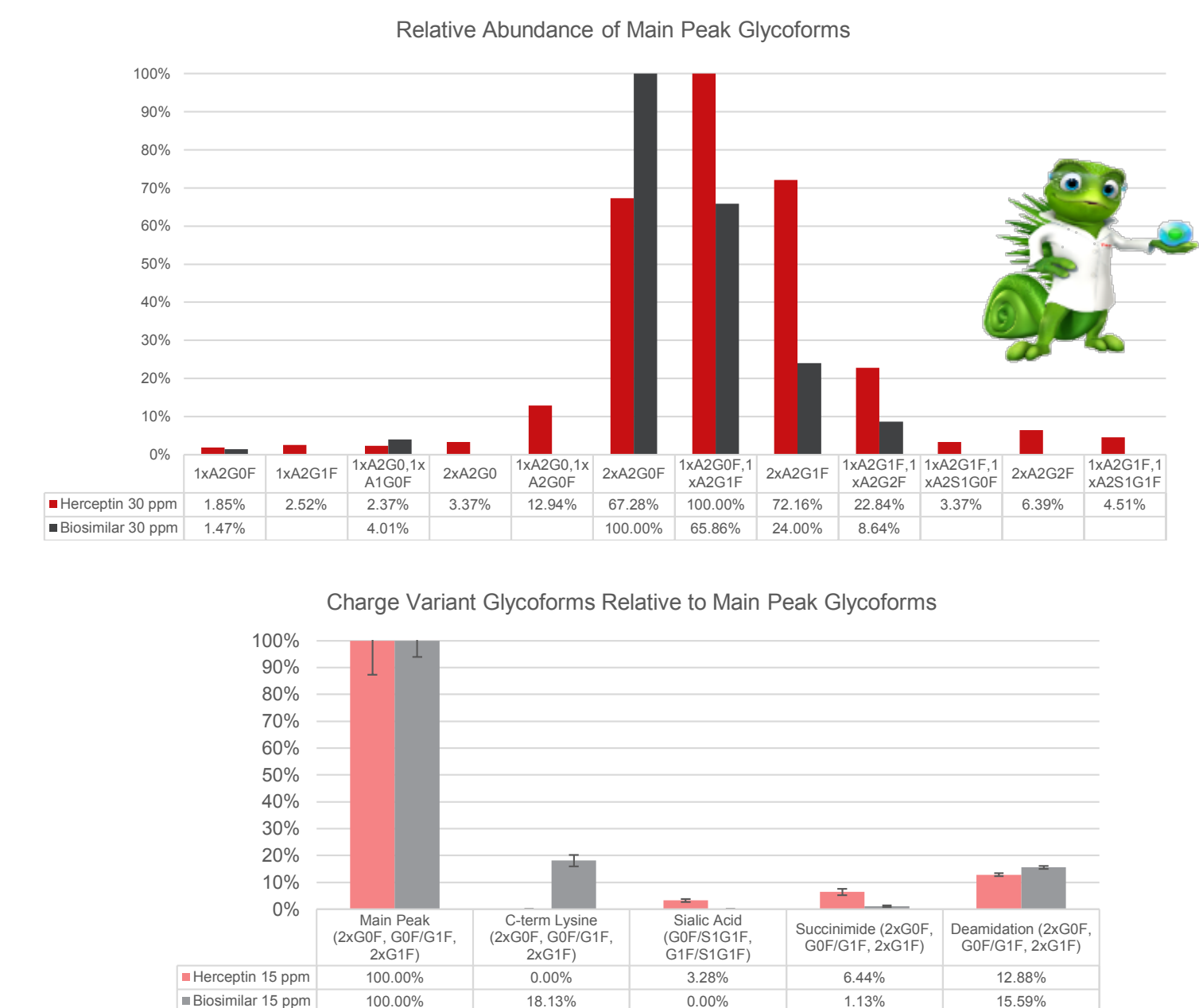


Figure 8. Intact glycoforms and charge variant CQAs can be simultaneously monitored. Charge variants CQAs are analyzed as groups of glycoforms, which may be useful for comparing innovator and biosimilar drugs. As shown below, we observed the biosimilar trastuzumab to contain decreased levels of succinimide isoforms compared to Herceptin, but to have markedly higher levels of C-terminal lysine retention.



CONCLUSIONS

- Method for monitoring multiple CQAs of an intact biotherapeutic compound to allow direct comparison of innovator and biosimilar products.
- Orbitrap mass analyzer is capable of mass measurements with high (sub-Dalton) precision
- BioPharma Finder software can be used to manually export mass lists from native ion exchange MS experiments, and m/z values may be targeted and quantified in Chromeleon software.

REFERENCES

- Harris RJ, Kabakoff B, Macchi FD, Shen FJ, Kwong M, Andeya JD, Shire SJ, Bjork N, Totpal K, Chen AB. Identification of multiple sources of charge heterogeneity in a recombinant antibody. J Chromatogr B Biomed Sci Appl. 2001 Mar;10(752(2)):233-45
- Rogers RS, Nightlinger NS, Livingston B, Campbell P, Bailey R, Ballard A. Development of a quantitative mass spectrometry multi-attribute method for characterization, quality control testing and disposition of biologics. MAbs. 2015;7(5):881-90. doi: 10.1080/19420862.2015.1069454.

TRADEMARKS/LICENSING

© 2018 Thermo Fisher Scientific Inc. All rights reserved. Herceptin is a registered trademark of Genentech, 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.

ThermoFisher
SCIENTIFIC