Characterization of Conformation of Therapeutic Antibody Aggregation with Optimized Hydrogen/Deuterium Exchange Mass Spectrometer

Terry Zhang¹, Shanhua Lin², Stephane Houel¹, Dave Horn¹, Xiaodong Liu², and Jonathan Josephs¹ ¹Thermo Fisher Scientific, San Jose, CA; ²Thermo Fisher Scientific, Sunnyvale, CA

ABSTRACT

Protein aggregation is one of the major critical quality attributes (CQA). To determine the structural change of antibody in connection with aggregation is important. In this study, a therapeutic antibody. Herceptin®, was treated with different acidic solutions (pH 0.5, 1.5 and 2.5) to induce aggregation. The aggregation was quantified by UV 280 absorption. The conformation of Herceptin and its aggregation was characterized by an optimized hydrogen/ deuterium exchange mass spectrometer (HDX). The HDX results revealed the more significant conformation change regions when aggregation was induced.

INTRODUCTION

Monoclonal antibodies (mAbs) have been increasingly used for detection and treatment of cancer and other diseases. Characterization of critical quality attributes (CQA) of mAb-based drugs is a primary concern for biopharmaceutical development. Structural characterization is used to assess the CQAs of biopharmaceutical products. Protein aggregation is a major CQA during monoclonal antibody (mAb) production. The formation of aggregates may impact safety and efficacy of mAbs. It is thus important to understand the mechanism of aggregation and the conformational changes of the aggregates. Hydrogen/deuterium exchange mass spectrometry (HDX) has emerged as a powerful tool to investigate the conformation of intact proteins, including mAbs. In this study, an optimized HDX workflow was used to probe the conformation of Herceptin and its aggregation.

MATERIALS AND METHODS

Instrument and Materials

A fully automated HDX platform, based on the H/D-X PAL[™] system (LEAP Technologies) and the Thermo Scientific[™] UltiMate 3000 pump system coupled online with an Thermo Scientific[™] UltiMate 3000 pump system coupled online with an Thermo Scientific[™] torbitrap Fusion[™] Thiofi[™] mass spectrometer, was used. Figure 1 shows the HDX work station. It is composed of an autosampler equipped with individually temperature-controlled sample plate and labeling plate. The syringe head and pepsin column can also be independently temperature-controlled. Chronos[™] control Software is fully integrated with the Thermo Scientific[™] Xcalibur[™] platform and the user interface is shown in Figure 2. Figure 3a) describes the three valve configurations at various stages of the experiment inside the cooling chamber. The flexibility allows for maximized productivity. Figure 3 b) shows the dual heads high pressure mixing pump on the left and ternary loading pump on the right. Flow rates and solvents can be changed on the fly during the experiment. The execution of the whole experiment the scentified and controlled by Chronos[™] Software.

Therapeutic antibody (Herceptin) aggregation was induced by adjusting the pH to 0.5, 1.5 and 2.5 with HCl and by incubating each sample at room temperature for 30 min. The pH of the samples was subsequently adjusted to pH 9. Both non aggregated and aggregated mAbs were diluted with labeling buffer and incubated for multiple time points. The samples were then quenched and digested online with a pepsin column in a fully automated manner using the H/D-X PAL system. The digested peptides were injected into a Thermo Scientific™ Acclaim™ Pepmap™ C18 reverse phase column with a 7 min gradient. MS analysis was performed with the Orbitrap Fusion mass spectrometer.

Data Analysis

Peptide identification was performed with Thermo Scientific™ Proteome Discoverer™ 1.4 software . Peptide mapping and PTM analysis was performed with Thermo Scientific ™ BioPharma Finder™ 1.0 software. HDX experimental data were analyzed with HDExaminer™ software (Sierra Analytics) and the HDX tool io ImAas Analyzer (licensed from Amgen).^{1,2}.





RESULTS

Herceptin and Herceptin Aggregation

Herceptin and its pH stressed samples were quantified by measuring UV 280 absorption. Figure 4 is the overlay absorption spectra of four samples. Aggregations around 25% for pH 0.5, 6% for pH 1.5 and 0% for pH 2.5 were observed.

Figure 2. Chronos software interface, integrated with Xcalibur software



Figure 3. a). Flexible three valve configuration for carrying out different experimental tasks simultaneously in the cooling chamber, b). NCS-3500RS with binary rapid separation micro flow pump with ternary loading pump



b

Figure 4. UV absorption spectrum of Herceptin and pH stressed samples



Figure 5. Peptide map of Herceptin by BioPharma Finder a) Heavy chain b) Light Chain





Peptide mapping of Herceptin

MS/MS experiments were first performed using non-deuterated samples for peptide identification. Nearly 100% sequence coverage was achieved for both the Herceptin and Herceptin pH stressed samples. Figure 5 is the peptide map of Herceptin generated by BioPharma Finder software. After online pepsin digestion, more than 200 and 100 peptides were respectively identified for heavy chain and light chains. These identified peptides were subsequently used to probe the conformation of the analyzed samples by HDX. Peptides modification summary was also generated by BioPharma Finder. For oxidation and deamidation, ther is no major difference between the control and pH stressed samples.



Herceptin and its Aggregation conformation characterization

Multiple time point HDX experiments were performed for both Herceptin and pH stressed Herceptin samples. Highly reproducible chromatograms were obtained for the various experimental time points. MS full scan spectra were collected to measure the deuterium uptake to probe the conformation.

Overall, the control and pH stressed samples had similar deuterium uptake profile but for most of the pH stressed sample peptides showed slightly more deuterium uptake, around 2 to 5%, than for the non-stressed sample. The differences could reach up to 10 to 15% for different regions for the light chain and could be even higher for the heavy chain. The sample stressed at pH = 0.5 showed more deuterium uptake than the pH = 1.5 sample. Finally, there was no significant deuterium uptake change between Herceptin and the stressed sample pH=1.5 (data not shown).

Figure 6. a): Herceptin and pH 0.5 stressed sample light chain deuterium uptake light chain residual plot. b), c): Example of specific peptides deuterium uptake plots of Herceptin and its pH 0.5 stressed sample



Figure 6 a) is the light chain deuterium uptake residual plot whereas figure b) and , c) are peptides uptake plots. In figure 6 b) peptide showed more differences for short labeling time but would reach to similar levels at longer time points, while in figure 6 c) peptide showed the significant differences persistent along all the time points. Protection factors were calculated and shown in figure 7. Certain regions have similar protection factors while other regions have significant differences, i.e., between amino acids160 and 190 for the heavy chain. These results were consistent with HDExaminer's analysis.

Figure 7. Herceptin and pH 0.5 stressed sample light chain protection factor comparison



Figure 8. Herceptin and pH 0.5 stressed sample heavy chain deuterium uptake mirror plots



Similar deuterium uptake profiles of Herceptin and its pH stressed sample were observed for heavy chain and are shown in figure 8. In general, the stressed sample showed more deuterium uptake, and the amount of uptake increased with degree of aggregation. This implies that aggregation would impact the 3D structure of Herceptin and more exposed to solvents, i.e., more deuterium uptake. Furthermore, the increase in deuterium uptake was not uniform across all the peptides. Example deuterium uptake dynamics and levels. With mass spectrometry, the regions with such increase can be pinpointed. A pronounced deuterium uptake increase (up to 40%, see left panel of figure 10) was observed from amino acid 120 to 200 and 240 to 255 of the heavy chain for the sample stressed at pH=0.5. The differences could be up to 5 daltons, see figure 10, right panel. These results indicated that aggregation would cause more conformational change for the Fab region

Figure 9. Herceptin and pH 0.5 stressed sample heavy chain peptides deuterium uptake plots







Figure 11. Herceptin and pH 0.5 stressed sample heavy chain protection factor comparison. Top: residual 1-200 Bottom: residual 201-449



To further understand the conformational dynamics, single amino acid residue level protection factors were calculated using Mass Analyzer HDX, (see figure 11). In most of the regions Herceptin showed slightly higher protection factor than the pH 0.5 stressed sample. From amino acid 110 to 210 and 240 to 255, Herceptin had a much higher protection factor, which is consistent with the deuterium uptake residual plots obtained from HDExaminer.

Figure 12. Herceptin and pH 0.5 stressed sample 1hr deuterium labeling Fab differential exchange map on crystal structure (PDB: 1N8Z)



Antibody aggregation is a complex, multistep process³. The conformation change with more deuterium uptake observed for the sample stressed at pH 0.5 agreed with the proposed aggregation mechanism^{3,4} and the reported results⁵ in the literature. The differential deuterium uptake of Herceptin and the sample stressed at pH 0.5 for1 hour labeling was mapped with Herceptin Fab crystal structure (PDB:1N82) and shown in figure 12. Most of the significant change regions were in the C_L, C_H and part of FC C_H2 domains.

CONCLUSIONS

 A fully automated HDX workflow was successfully applied to the study of conformational changes of Herceptin upon aggregation.

-In this study, we were able to pinpoint the subtle but significant changes around the $C_{\rm L},\,C_{\rm H}$ and $C_{\rm H}2$ regions.

•The MS data were analyzed by two independent software packages (HDExaminer and Mass Analyzer) and the conclusions were consistent.

REFERENCES

- 1. 1. Mass Analyzer, Amgen Inc.
- 2. Z. Zhang et al, Anal. Chem. 2012, Vol. 84, 4942-493
- 3. D. Weis et al, Hydrogen Exchange Mass Spectrometry of Proteins
- 4. P. Arosio et al, Pharmaceutical Research, 2013, Vol. 30, 641-654
- 5. R.E. lacob, et al, J. Pharmaceutical Science, 102,(12), 4315-4329

www.thermofisher.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. Herceptin is a trademark of Genetech Inc. HD/X-PAL and Chronos are trademarks of LEAP Technologies. HDExaminer is a trademark of SierraAnalytics. ISO is a trademark of the International Standards Organization. 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 233555 USA +1 800 532 4752 PN64771-EN 0616S



A Thermo Fisher Scientific Brand