

Imaged capillary isoelectric focusing – Mass Spectrometry (iCIEF- MS) online coupling for polatuzumab vedotin charge heterogeneity analysis using native MS

Xiaoxi Zhang¹, Tony Chen², Tao Bo², Tiemin Huang², Min Du³.

¹ Thermo Fisher Scientific, Shanghai, China, ² Advanced Electrophoresis Solutions Ltd., Cambridge, Canada, ³ Thermo Fisher Scientific, Lexington, MA, US

Abstract

Purpose: To demonstrate the capability of online coupling iCIEF-MS for cysteine-linked ADC charge variant analysis under near native condition.

Methods: CEInfinite iCIEF platform (AES Ltd., Canada) and Thermo Scientific™ Orbitrap Exploris™ 240 mass spectrometer with BioPharma option were employed.

Results: Successfully separated and identified charge variants of polatuzumab vedotin, a latest-generation cysteine-linked ADC using iCIEF-MS online coupling platform under native condition.

Introduction

Antibody-drug conjugates (ADCs) represent the forefront of the next generation of biopharmaceuticals. An ADC typically comprises an antibody covalently linked to a cytotoxic drug via a linker, resulting in a highly heterogeneous product. The conjugation will bring extra charge heterogeneity to the biomolecule, therefore, charge based separation technologies are important in ADC characterization.

iCIEF has become an indispensable tool in therapeutic protein development and manufacturing because of its high analytical throughput, ease of use, fast method development and excellent reproducibility. Recently, iCIEF-MS has attracted much attention to utilize for protein charge variant analysis. Here we did online iCIEF-MS to analyze charge heterogeneity of polatuzumab vedotin (Polivy®, Roche), a latest-generation cysteine-linked ADC, using CEInfinite iCIEF platform and Thermo Scientific™ Orbitrap Exploris™ 240 mass spectrometer. NISTmAb was used for method development for iCIEF-MS online coupling under near native condition.

Materials and methods

Sample Preparation

Commercially available polatuzumab vedotin was dissolved in ddH₂O and desalted using 10k cut-off filter. Final protein concentration is 1.5mg/mL in ddH₂O with 1% HR3-10 and 1% HR6-8 carrier ampholytes.

iCIEF Focusing and mobilisation

200 µm ID acrylamide derivative coated (AD) capillary cartridge was used. The focusing was performed using 1 min-1000 V, 1 min-2000 V and 10 min-3000 V. 3000 V was applied during mobilisation of focused protein bands; the mobilisation speed was 70 nL/min with 0.1% FA, H₂O and 3 µL/min make up solution (10mM NH₄Ac:ACN=9:1) added through a micro tee. Mobilisation time was 50 min.

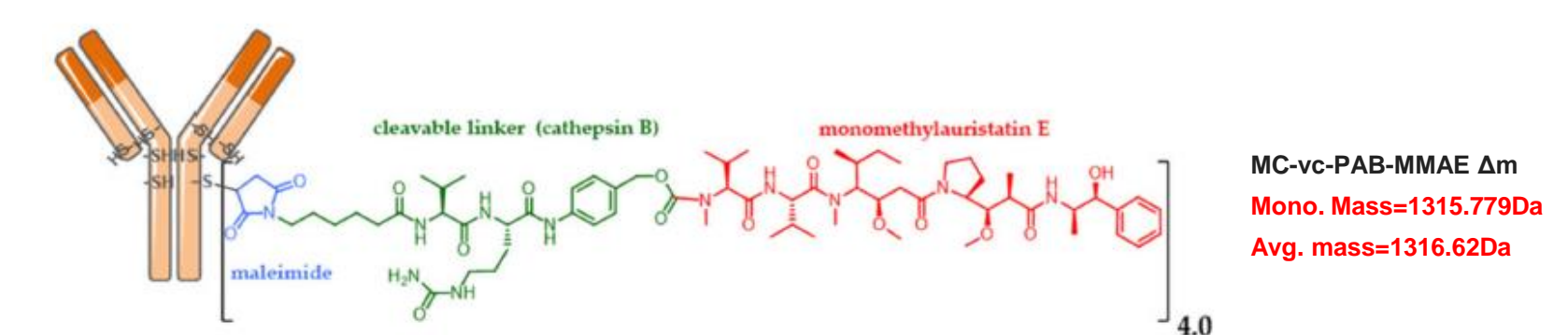
Mass Spectrometer:

An Orbitrap Exploris™ 240 mass spectrometer with BioPharma option was used for data acquisition. Both iCIEF and MS platforms were controlled using Thermo Scientific™ Chromeleon™ software. MS settings are listed in Table 1.

Data Analysis

Data analysis was performed using Thermo Scientific™ BioPharma Finder™ software 5.2.

Figure 1. The Schematic of polatuzumab vedotin.



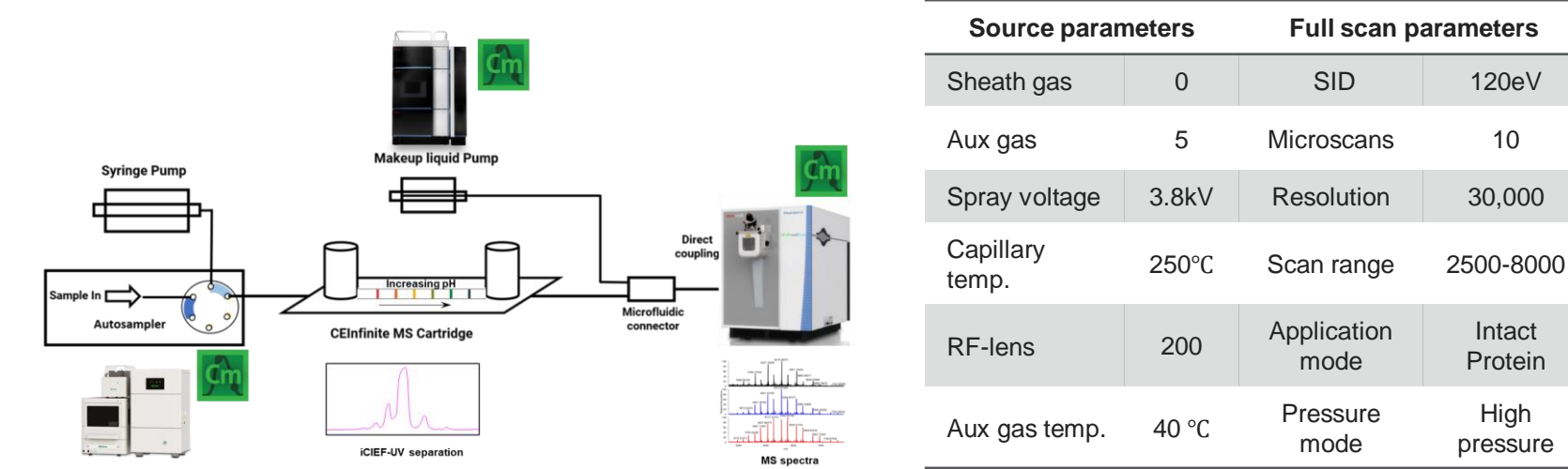
Results

Method development and optimization for iCIEF-MS online coupling under near native condition

Polatuzumab vedotin, a latest-generation cysteine-linked ADC was used in this study. The schematic of polatuzumab vedotin is shown in Figure 1^[1]. During the conjugation of polatuzumab vedotin, the interchain disulfide bridges of mAb were partially reduced, and the mAb was conjugated to a maleimidocaproyl-valinecitruilline-p-aminobenzyloxycarbonyl linker- monomethyl auristatin E (MC-vc-PAB-MMAE), leads to a 0/2/4/6/8 drug payload distribution. The average DAR of this ADC is around 4 according to previous publication^[1]. As the interchain disulfide bridges of the mAb were reduced for conjugation, the light and heavy chains (LC and HCs) are non-covalently bonded. Therefore, native intact MS is essential for charge heterogeneity analysis of polatuzumab vedotin at the intact ADC level.

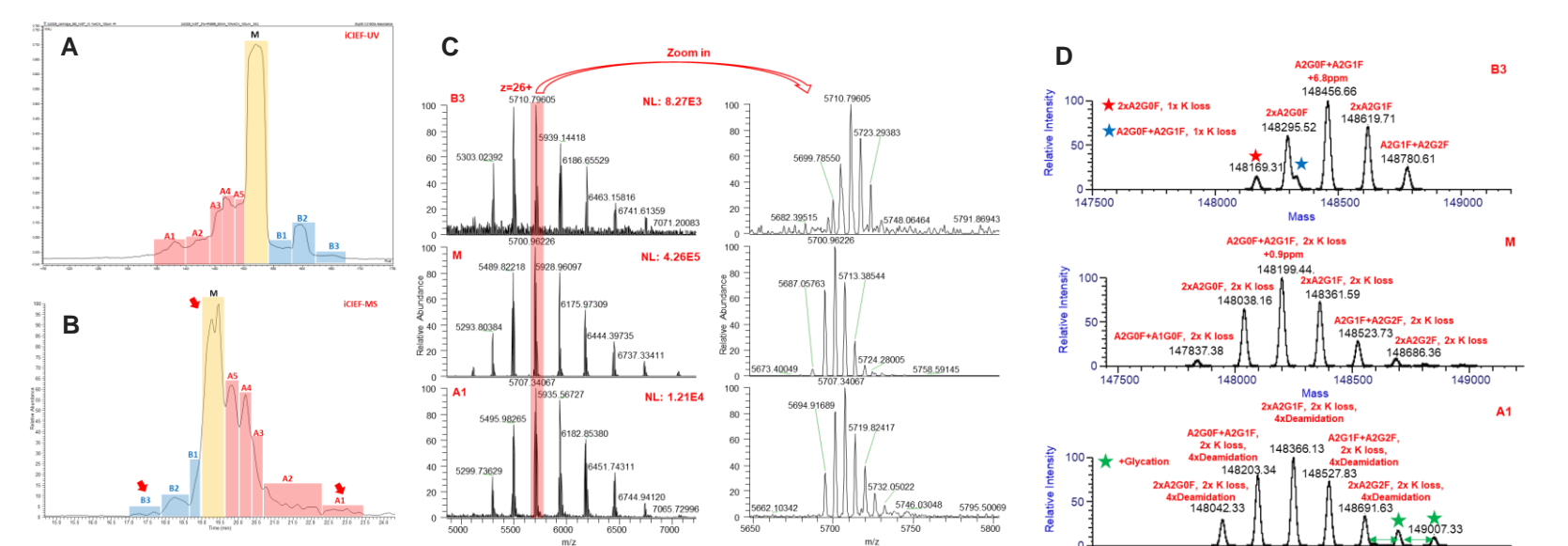
Figure 2 illustrates iCIEF-MS online coupling workflow. After proteins are separated into groups based on their isoelectric point (pI), the subsequent mobilization process introduces protein bands into MS to isolate the charge variant peaks. During pressure mobilization, an electric field keeps the samples in the separation capillary focused. Thermo Scientific™ Chromeleon™ software can control CEInfinite iCIEF platform and Thermo Scientific™ Orbitrap Exploris™ 240 mass spectrometer for data acquisition, which better meets compliance requirements.

Figure 2. iCIEF-MS online coupling workflow.



iCIEF-MS online coupling under denaturing condition is a mature workflow now and has been used in mAb and fusion protein charge variant analysis^[2-4]. However, iCIEF-MS online coupling under native condition is quite challenging because the MS signal intensity of protein is much lower under native condition compares to denaturing condition, which means the sensitivity of MS is critical, especially for low abundance charge variants identification. Also, the make-up liquid needs to be changed to MS compatible buffer while keeping samples stay native. Figure 3 shows the iCIEF-MS online coupling under native condition of NISTmAb charge variant analysis. It is clearly that charge variant peaks were well separated by iCIEF, benefited from highly sensitivity of Orbitrap analyzer, even the low abundance basic peak (B3) components had high S/N in raw MS spectra. The mass error of A2G0F+A2G1F, 2x K loss in main peak is 0.9ppm, demonstrating the highly mass accuracy of Orbitrap platform.

Figure 3. The charge variant analysis of NISTmAb (1.6µg sample was loaded) using online coupling iCIEF-MS under near native condition. A, iCIEF-UV profile. B, iCIEF-MS profile. C, raw spectra of basic peak3(B3), main peak(M) and acidic peak1(A1), as labelled in 3B. D, deconvoluted spectra of B3, M and A1.



iCIEF-MS online coupling of polatuzumab vedotin under near native condition

As discussed in previous section, to keep the protein in native state, the make-up liquid and mobilisation liquid composition are critical. For native mAb analysis, the combination of make-up liquid (10mM NH₄Ac:ACN=9:1) and mobilisation liquid (0.1%FA, H₂O) can maintain mAb at near native condition, as shown in Figure 3. However, when preparing ADC sample for iCIEF analysis, a common approach is to add formamide into sample solution to improve solubility. But formamide will break the non-covalently bonds between chains of cysteine-linked ADC. Figure 4 displays iCIEF-UV profile and MS spectra of main peak, with 0%, 1% and 5% formamide in polatuzumab vedotin (v/v) respectively. It is clearly that formamide leads to better peak shape and separation at iCIEF level, but even with lower to 1% formamide, the LC/HC dissociation can be observed in MS spectra. To keep the ADC sample at native state, we choose no formamide in ADC sample as the final condition. iCIEF-MS online coupling of polatuzumab vedotin was performed with optimized experiment make-up liquid (10mM NH₄Ac:ACN=9:1), mobilisation liquid (0.1%FA, H₂O) and no formamide in sample. Figure 5 A and B display iCIEF-UV and iCIEF-MS profiles, four acidic peaks, main peak and two basic peaks were separated and identified.

Figure 4. Sample buffer optimization for polatuzumab vedotin.

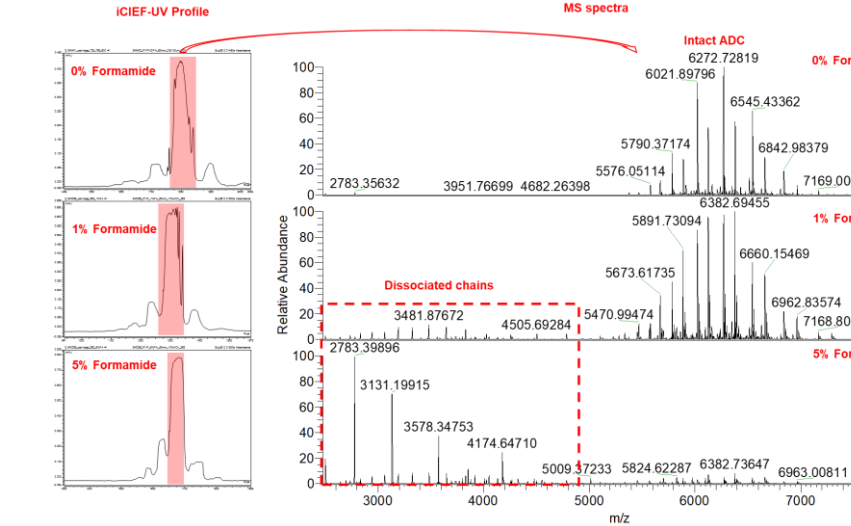
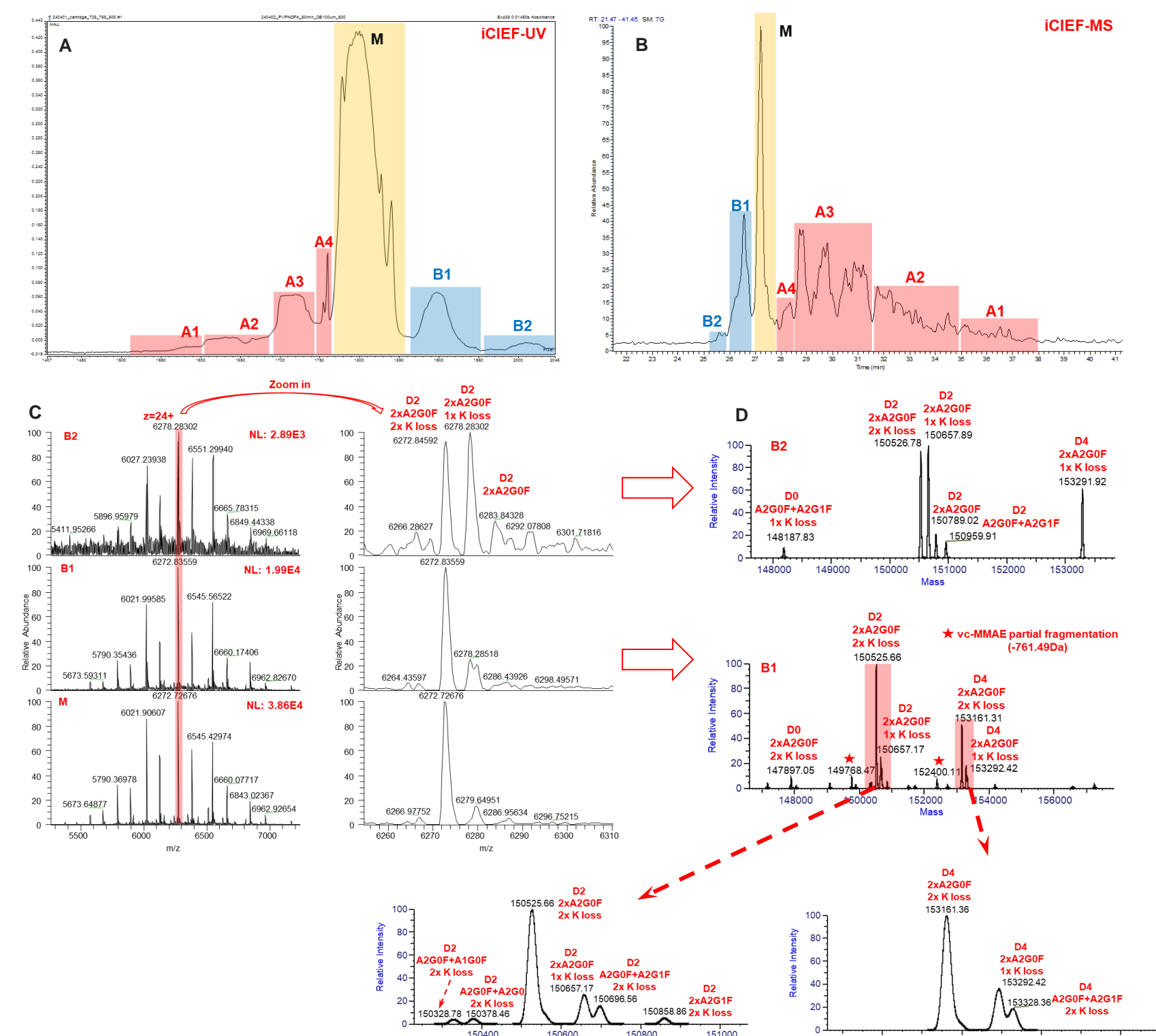
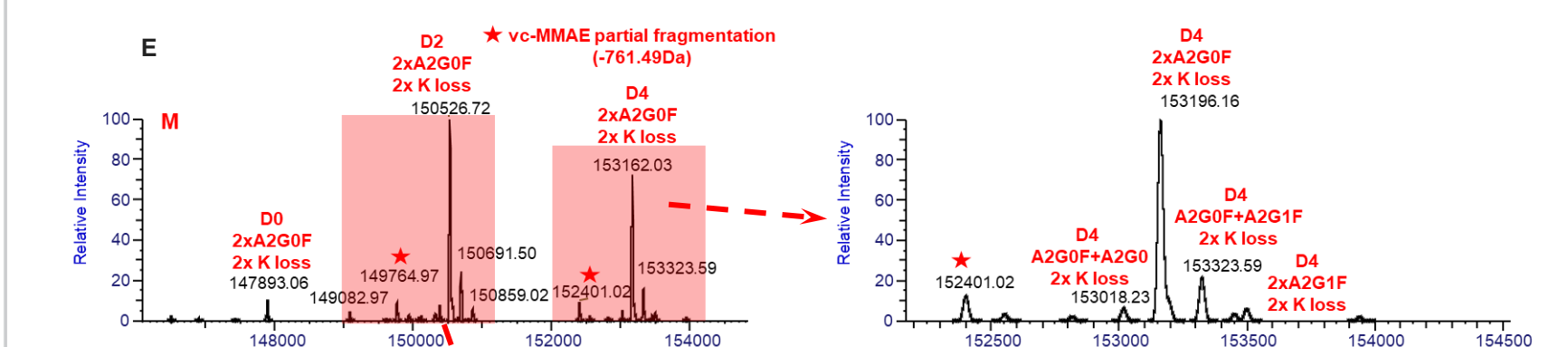


Figure 5. The charge variant analysis of polatuzumab vedotin (2.4µg sample was loaded) using online coupling iCIEF-MS under near native condition. A, iCIEF-UV profile. B, iCIEF-MS profile. C, raw spectra of basic peak1-2 (B1-2) and main peak (M). D, deconvoluted spectra of B1 and B2. E, deconvoluted spectra of main peak. F, raw spectra of acidic peak1-4 (A1-4). G, deconvoluted spectra of acidic peak1-4 (A1-4).

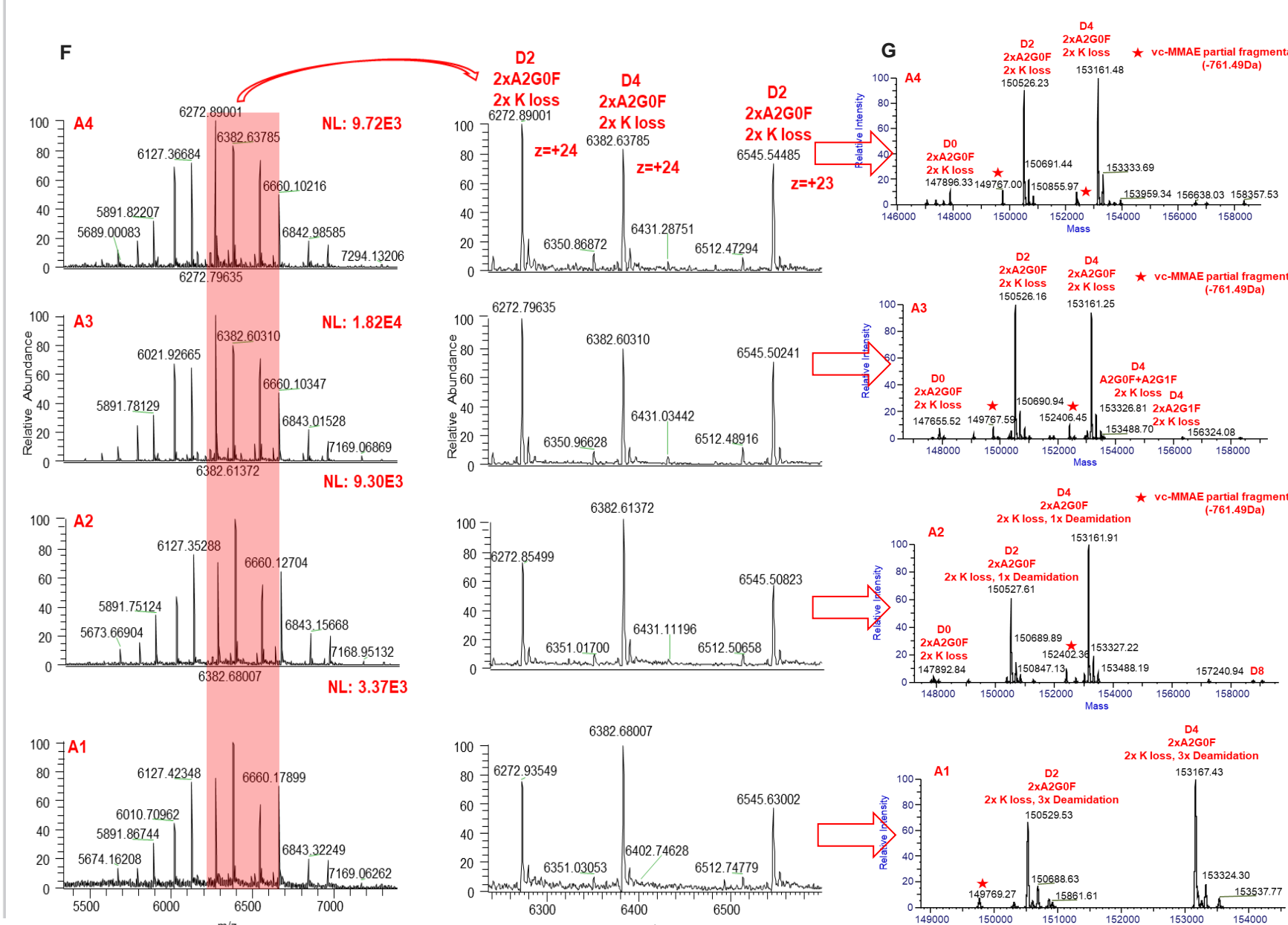


Without formamide in sample, we lost separation resolution at iCIEF level (Figure 5A), but Orbitrap based MS platform could provide high resolution to compensate the loss. Another advantage of Orbitrap MS platform is excellent sensitivity. In this experiment, only 2.4µg sample was loaded, and compares to denaturing condition, the MS signal of intact protein is much lower under native condition. Even for low abundance basic peak 2 (relative abundance 0.79%), the S/N of raw spectra is high (Figure 5C). D2 and D4 carrying one or two lysine at C-terminal of heavy chain were identified in B2 (Figure 5D). B1 is mixture of D2 and D4 with 2x K loss or 1x K loss. The N-glycosylation distribution of D2 and D4 were shown in expanded view of Figure 5D. Low abundance N-glycoforms, such as A2G0F+A1G0F and A2G0F+A2G0 were detected. Vc-MMAE partial fragmentation mass shift (-761.49Da) was observed and labelled with red star.



The major components in main peak are D2 and D4 N-glycosylation distribution with 2x K loss (Figure 5E). Figure 5F displays the raw spectra of acidic peaks. In acidic peak A4 and A3, D2/D4 ratio varied while components are almost the same; the deamidated variants were detected in acidic peak A2 and A1. With increasing pI, the relative abundance% of D4 increased. In basic and main peak, the top drug payload isoform is D2 while in acidic peak A4, A2 and A1, D4 became the top drug payload isoform, indicates drug conjugation may affects charge heterogeneity of ADC, especially the net charge distribution. It looks like iCIEF separates charge variants of cysteine-linked ADC based on PTM-induced charge heterogeneity more than bioconjugation-induced heterogeneity.

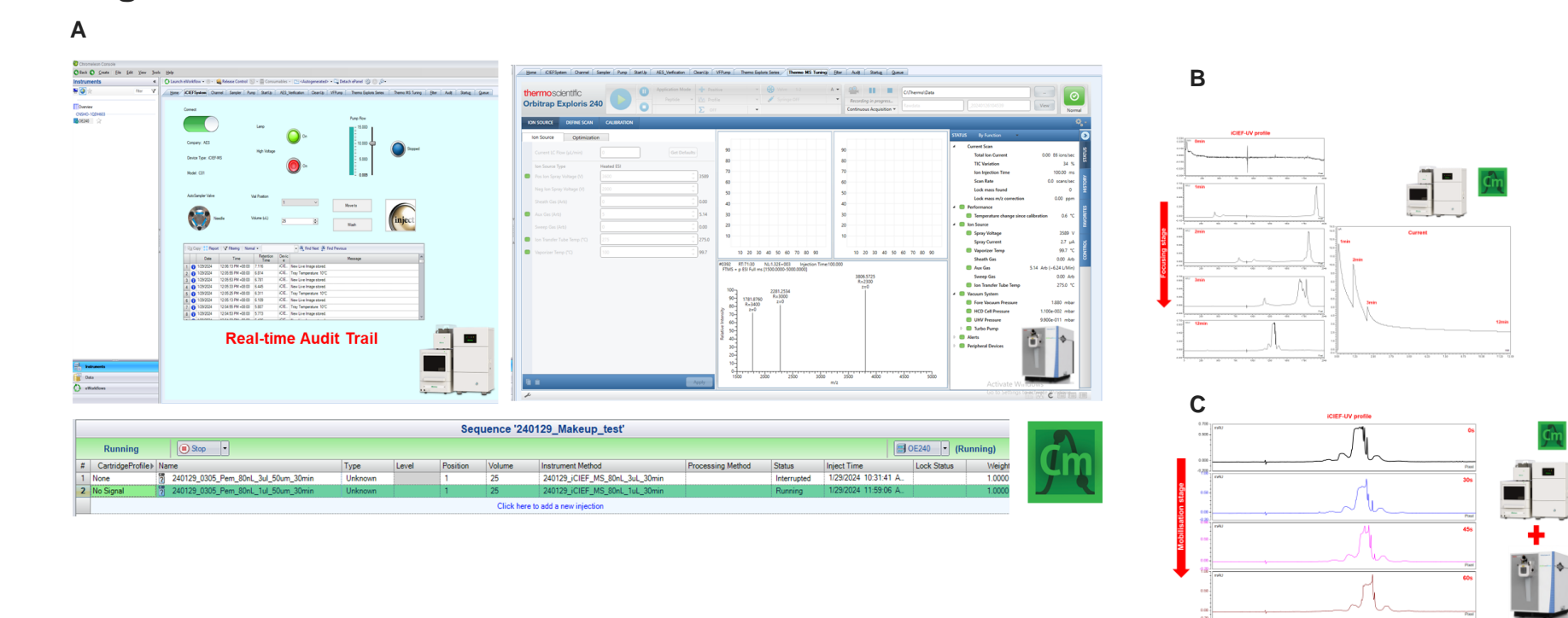
The major components in main peak are D2 and D4 N-glycosylation distribution with 2x K loss (Figure 5E). Figure 5F displays the raw spectra of acidic peaks. In acidic peak A4 and A3, D2/D4 ratio varied while components are almost the same; the deamidated variants were detected in acidic peak A2 and A1. With increasing pI, the relative abundance% of D4 increased. In basic and main peak, the top drug payload isoform is D2 while in acidic peak A4, A2 and A1, D4 became the top drug payload isoform, indicates drug conjugation may affects charge heterogeneity of ADC, especially the net charge distribution. It looks like iCIEF separates charge variants of cysteine-linked ADC based on PTM-induced charge heterogeneity more than bioconjugation-induced heterogeneity.



Chromeleon control for both CEInfinite iCIEF platform and Orbitrap Exploris™ 240 mass spectrometry

Since the driver of CEInfinite iCIEF platform has been integrated into Chromeleon from 7.3.1, Chromeleon can now be used for instrument control of CEInfinite coupled to multiple Orbitrap series MS. Figure 6A shows Chromeleon software page, including CEInfinite iCIEF platform page, Thermo MS tuning page and running sequence. Both real-time and historical audit trails log can be read or exported according to users needs. Figure 6B is iCIEF-UV profiles and current channels during focusing stage and Figure 6C is iCIEF-UV profiles during mobilisation stage. All information were recorded in Chromeleon data file to meet the regulatory requirements.

Figure 6. Chromeleon control for both CEInfinite iCIEF platform and Orbitrap Exploris™ 240 mass spectrometer. A, software interface. B, iCIEF-UV profiles and current channels during focusing stage. C, iCIEF-UV profiles during mobilisation stage.



Conclusions

- In this work, we developed and optimized online coupling iCIEF-MS methods for charge variant analysis of NIST mAb and polatuzumab vedotin under near native condition.
- Using NISTmAb, we demonstrated the developed method was highly sensitive and provide excellent mass accuracy.
- We observed the iCIEF based charge variants profile of polatuzumab vedotin was more influenced by the mAb charge heterogeneity (such as K variants and PTMs) than conjugation-induced heterogeneity.
- In addition, we successfully demonstrated the feasibility of Chromeleon control for CEInfinite iCIEF platform coupled to Orbitrap Exploris™ 240 mass spectrometry.

References

- Pharmaceuticals 2020, 13, 245; doi:10.3390/ph13090245.
- Analytica Chimica Acta, 1297 (2024) 342349. <https://doi.org/10.1016/j.aca.2024.342349>.
- Analytical Biochemistry, Volume 680, 1 November 2023, 115312;
- Analytical Biochemistry, Volume 660, 1 January 2023, 114961.

Trademarks/licensing

© 2024 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. PO253-2024-EN