

# Capillary Electrophoresis – Mass Spectrometry for Intact Mass Analysis of Antibodies and Antibody-Drug-Conjugates

Michael Wenz<sup>1</sup>, Andreas Krupke<sup>1</sup>, Chien-Hsun Chen<sup>2</sup>, Stephane Houel<sup>2</sup>, Daniel Lopez-Ferrer<sup>2</sup> and Aran Paulus<sup>2</sup>

<sup>1</sup>Thermo Fisher Scientific, Life Sciences Solutions, South San Francisco, CA 94080; <sup>2</sup>Thermo Fisher Scientific, Chromatography and Mass Spectrometry, San Jose, CA 95134

## ABSTRACT

Correct identification of heterogeneity of monoclonal antibodies (mAbs) and antibody-drug-conjugates (ADCs) is an essential analytical tool in the biopharmaceutical industry. High resolution accurate mass spectrometry has established itself as a powerful tool for the characterization of such compounds; however some proteoforms might not be observed due to potential sample complexity. Here, we used the ZipChip system from 908 Devices made of micro-fabricated glass chips, coupled to a Thermo Scientific™ Q Exactive™ Biopharma system to separate and characterize charge variants of the NIST mAb standard and an ADC derived from it. We were able to identify the three published charge variants with 0, 1 or 2 lysine residues. Using the Thermo Scientific™ Biopharma Finder™ 2.0 software package, accurate intact mass for at least 15 glyco-forms were obtained. The NIST mAb has only one glycan site per heavy chain, which was modified using click chemistry to generate an ADC. The resulting model ADC also had charge variants, which were successfully separated and annotated with the correct MS1 by CEMS.

## INTRODUCTION

The main goal of this work was to develop a method to characterize heterogeneous mAb and ADCs in a top down approach. This was achieved by combining the powerful combination of efficient separation by capillary electrophoresis (CE) and HRAM together with data analysis software. The separation of proteoform charge variants is quite a challenge using chromatography, but CE provides an ideal separation technique for this problem. Microfluidic CE is advantageous for the mAb analysis given its good separation efficiency and short migration times because on high field strength and short separation channel length. The Orbitrap mass spectrometer family is also a good option for this application, and the Thermo Scientific™ Q Exactive™ HF mass spectrometer with BioPharma option and extended mass range of up to 8000 D is in particular suitable for the analysis of very large biomolecules. The BioPharma Finder software package is specifically designed to characterize complex proteins using dedicated deconvolution of mass spectra for individual time windows in combination with a thorough comparison to a reference database with all variants. In this study, a powerful combination of microfluidic CE separation, MS detection, and data analysis is presented to solve one of the most challenging workflows in the Biopharma industry.

## MATERIALS AND METHODS

The NIST reference material (NIST mAb standard RM 8671) used in this study has a published sequence and post translational modifications (PTMs). It is a recombinant humanized IgG1k antibody with any process-related impurities removed through various purification steps. The heavy chain of this mAb is known to have a high abundance of PTMs such as N-terminal pyroglutamination, c-terminal lysine clipping, and glycosylation. 10µg/µL of raw sample was diluted with deionized water to 0.5µg/µL prior to injection. No desalting procedure is necessary for this series of CEMS experiments. The HR ZipChip was primed with background electrolyte solution (BGE) consisting of 0.2% acetic acid and 10% IPA (pH=3.17). Once a stable electrospray was observed, 10µL of 0.5µg/µL mAb was pipetted into the sample reservoir. Pressure injection with 2 psi for 4s was used for sample loading, introducing 0.4nL corresponding to 0.2ng of mAb into the separation channel. High voltage of 20kV was applied for CE separation within 22 cm-length of channel, and ~2kV was used for nano-ESI. Each run was finished in less than 3 minutes.

The NIST mAb standard was labeled using click chemistry with a negatively charged fluorescein derivative or biotin, which were attached on both N-linked glyco units on the heavy chain, adding either two times 1145.37 D for AF488 or two times 857.03 D for the biotin label to the MW of the antibody. The Q Exactive HF Biopharma was set to the optimized conditions for ultra-high mass biomolecule, and to maintain a scan rate of 6 scans/sec. In-source CID (100 e.v.) was applied to assist desolvation. m/z range (2500–6000) was selected for this mild-denatured condition. BioPharma Finder 2.0 was used in this study for data analysis. The sequence and PTM information of the NIST mAb was first set up in protein sequence manager. The identification of different variants can be accomplished after the mass spectra were deconvoluted to the intact mass with ReSpect.



Figure 1. Characterization of mAb/ADC is accomplished by separation with the ZipChip device, on-line analysis with high resolution Q Exactive Biopharma mass spectrometer, and identification with BioPharma Finder software.

## Instrumentation

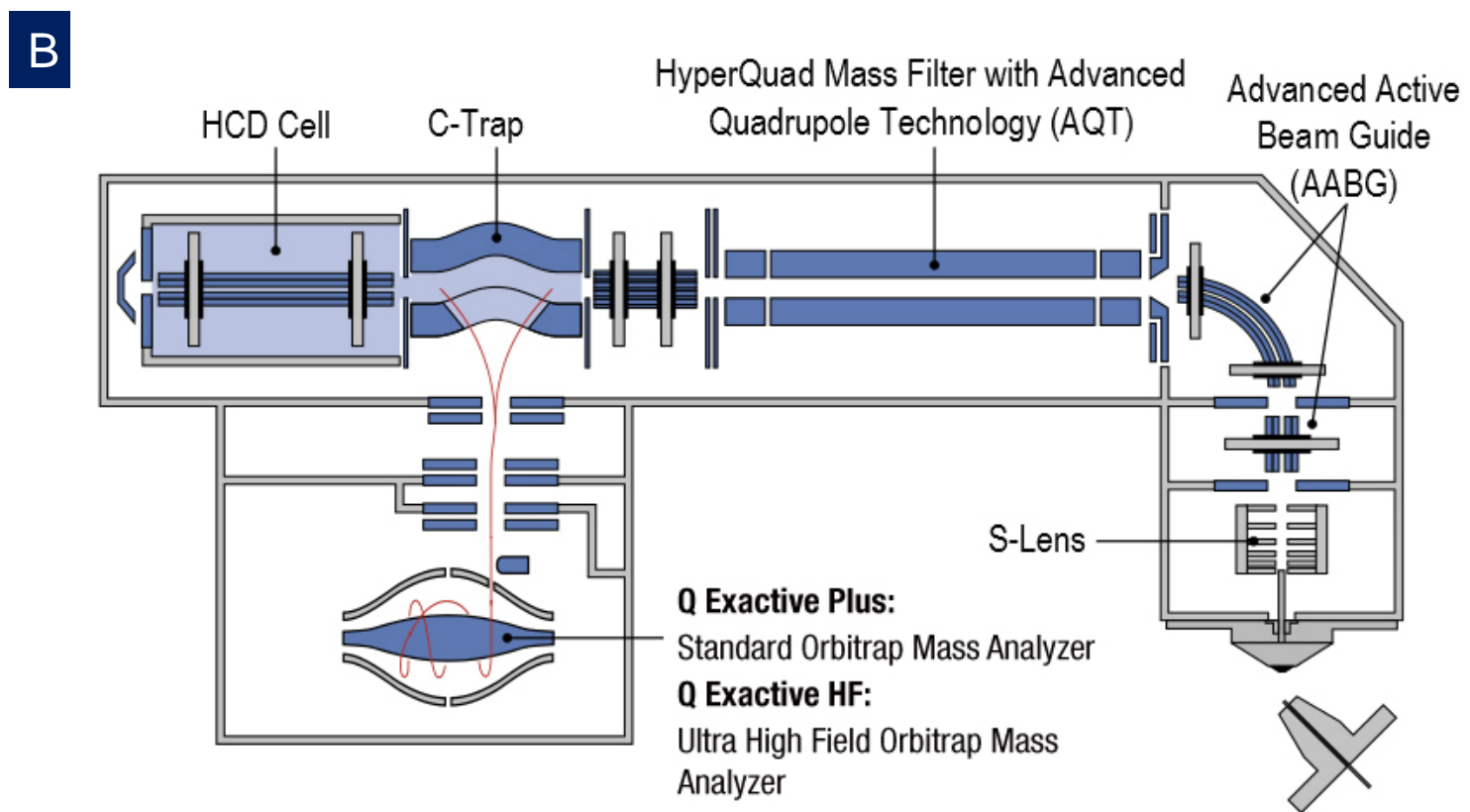
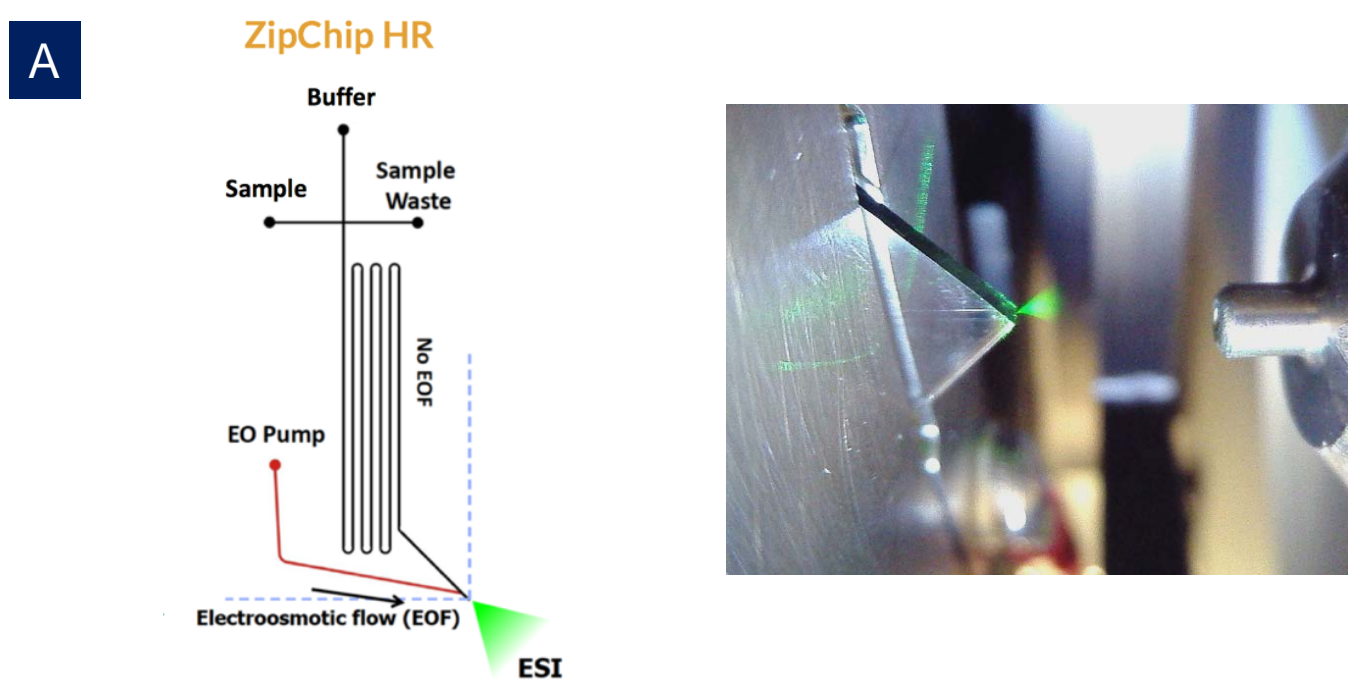


Figure 2. A) Schematic of high-resolution (HR) microfluidic CE chip with a 22 cm channel and integrated nano-ESI emitter. The nano-ESI spray can be monitored with the integrated green laser and camera. B) Schematic of Q Exactive HF mass spectrometer including the high-performance quadrupole precursor selection and high-resolution, accurate-mass orbitrap detection. The BioPharma Finder software package is optimized for 3 protein workflows: high-mass-range(HMR) mode for intact analysis, intact protein mode for subunit analysis, and normal mode for peptide mapping.

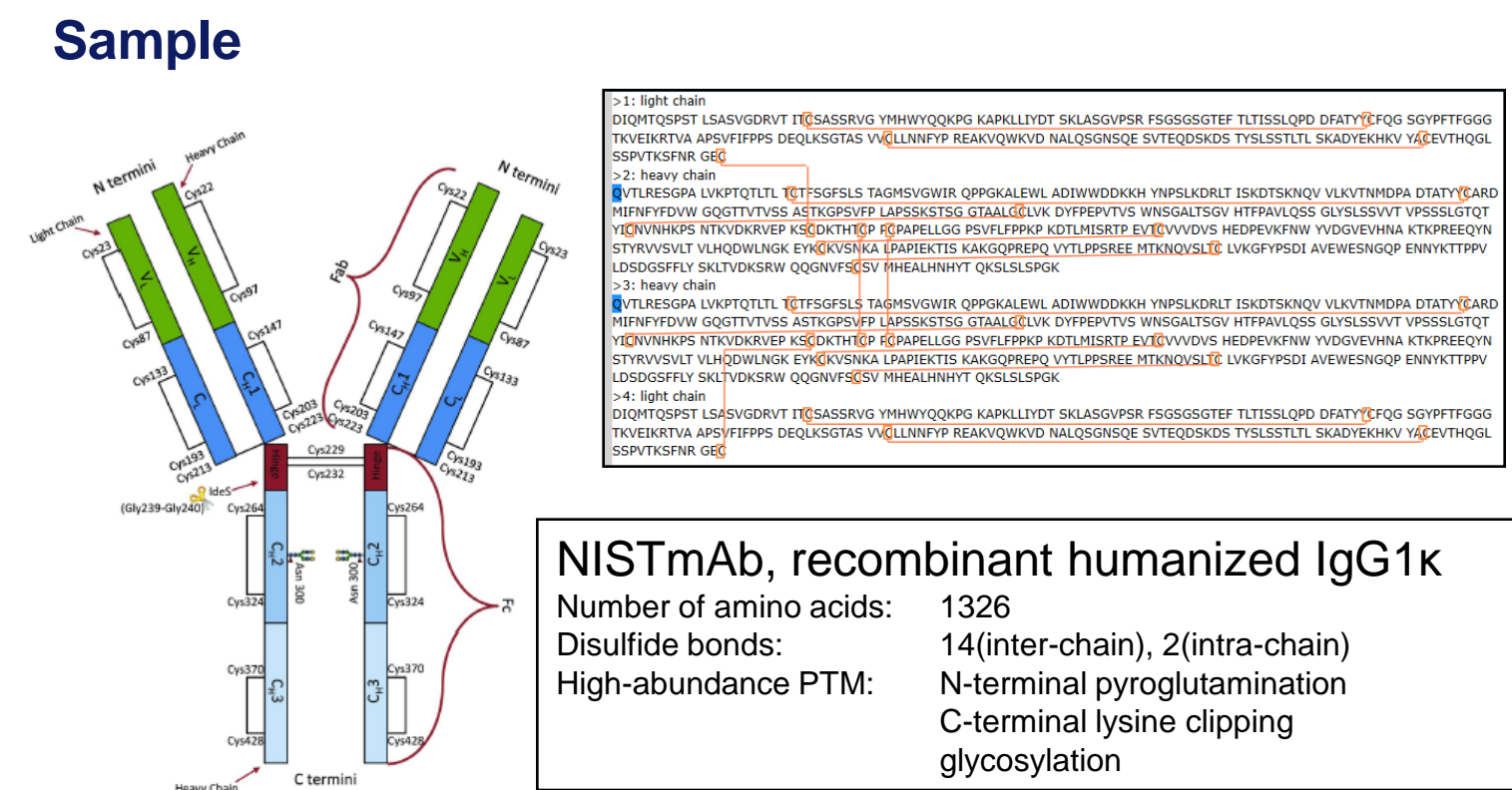


Figure 3. NISTmAb 8671 including amino-acid sequence, disulfide-bond linkage, and post-translational modification.

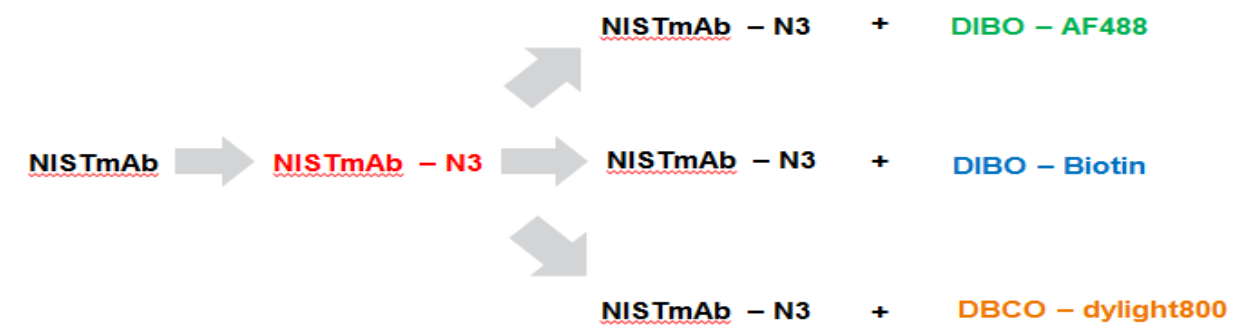


Figure 4. Reaction scheme to obtain ADCs from the NIST mAb standard using click chemistry. Three different payloads were generated, using AF488, Biotin and dylight800. CEMS analysis in Figure 8

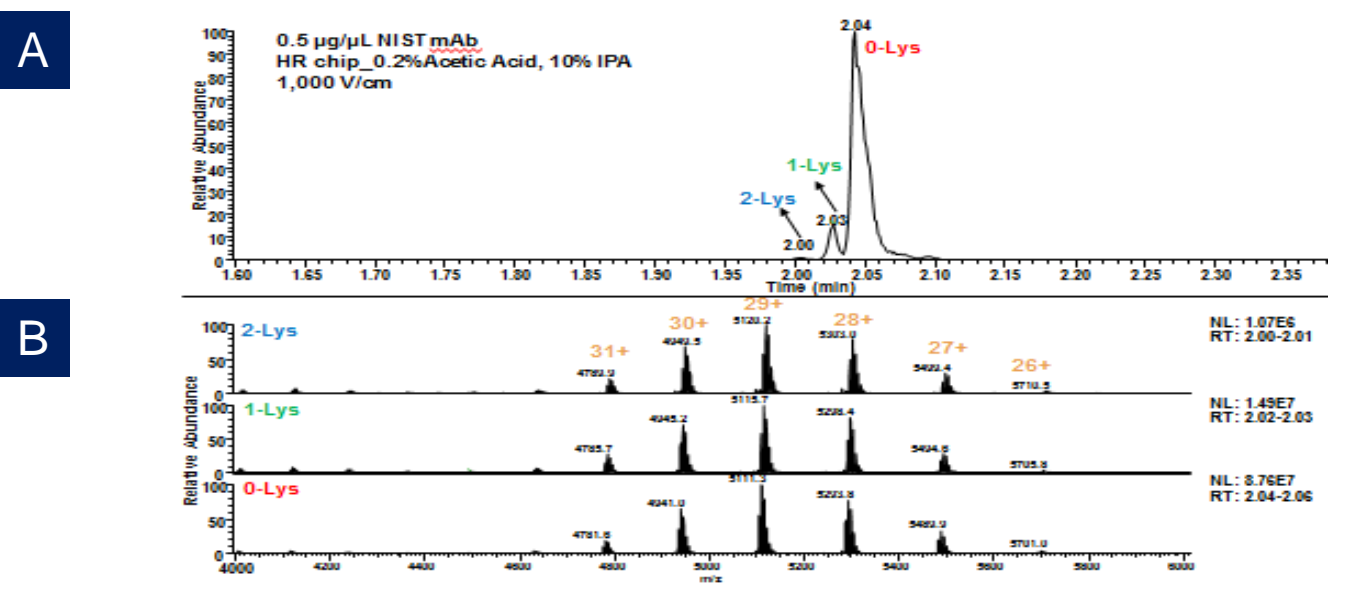


Figure 5. A) Electropherogram of the NIST mAb in a 2.35 min run on a HR microfluidic chip hyphenated to Q Exactive Biopharma mass spectrometer. Three different lysine variants were separated. B) Three mass spectra were extracted from the three peaks of the electropherogram. Lysine variant (128 Da) can be identified after the comparison of deconvoluted mass spectra. Higher number of lysine on the C-terminal of heavy chain provides higher mobility due to the additional positive charges from lysine.

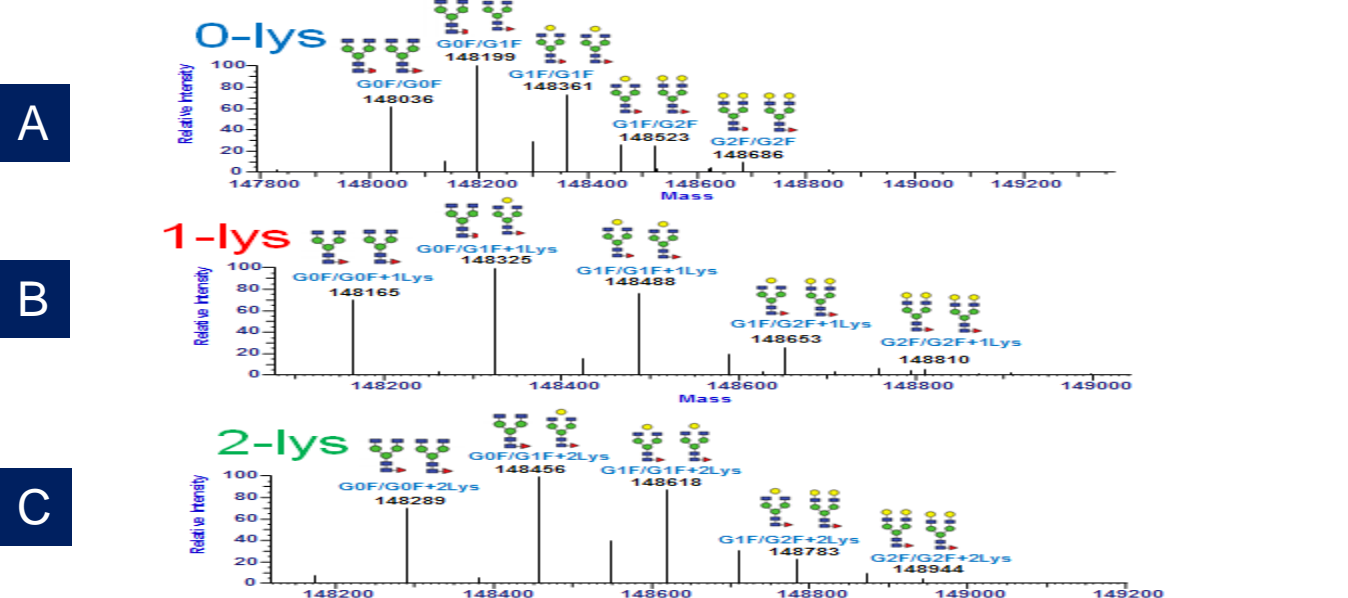


Figure 6. Three deconvoluted mass spectra from different migration time intervals, corresponding to different lysine variants. A) mAb with 0 lysine adduct including 5 glycosylated variants was detected B) mAb with 1 lysine adduct including 4 glycosylated variant was detected C) mAb with 2 lysine adducts including 5 glycosylated variants were detected.

Protein Name	Modification	Average Mass	Theoretical Mass	Mass Error (ppm)	Relative Abundance	RT Range
1 NIST mAb	1xG0F1F	148199.98	148199.21	6.3	100.00	2.036-2.078
2 NIST mAb	1xG1F0F1	149361.07	148361.38	1.9	99.33	2.036-2.078
3 NIST mAb	1xG0F0F	149036.73	148037.07	11.2	60.90	2.036-2.078
4 NIST mAb	1xG1F02F	148522.19	148523.48	0.7	44.32	2.036-2.078
5 NIST mAb	1xG2F02F	148852.99	148885.65	20.8	10.78	2.037-2.051
6 NIST mAb plus 1K	1xG0F1F	148327.28	148327.41	0.9	10.67	2.013-2.036
7 NIST mAb plus 1K	1xG1F0F1	148490.08	148488.55	3.6	8.32	2.013-2.036
8 NIST mAb plus 1K	1xG0F0F	148166.99	148166.27	4.9	6.65	2.021-2.034
9 NIST mAb plus 1K	1xG1F02F	148651.79	148651.68	0.7	4.28	2.017-2.036
10 NIST mAb plus 1K	1xG2F02F	148811.51	148813.83	16.5	0.80	2.017-2.026
11 NIST mAb plus 2K	1xG1F0F1	148619.07	148617.75	8.9	0.67	1.995-2.013
12 NIST mAb plus 2K	1xG0F1F	148456.13	148456.61	3.2	0.53	2.001-2.012
13 NIST mAb plus 2K	1xG0F0F	148294.46	148293.47	6.7	0.46	1.996-2.012
14 NIST mAb plus 2K	1xG1F02F	148789.80	148779.88	6.1	0.33	1.999-2.012
15 NIST mAb plus 2K	1xG2F02F	148942.99	148942.03	6.4	0.07	2.003-2.011

Figure 7. Identified proteoforms of the NIST mAb through the BioPharma Finder 2.0 software.

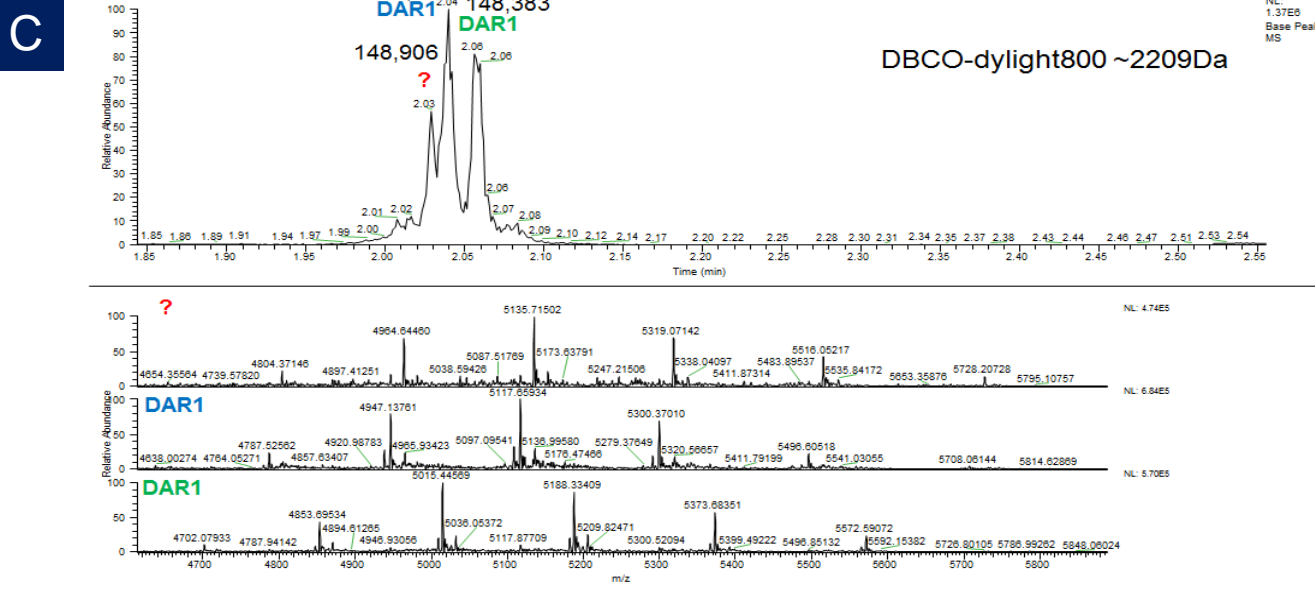
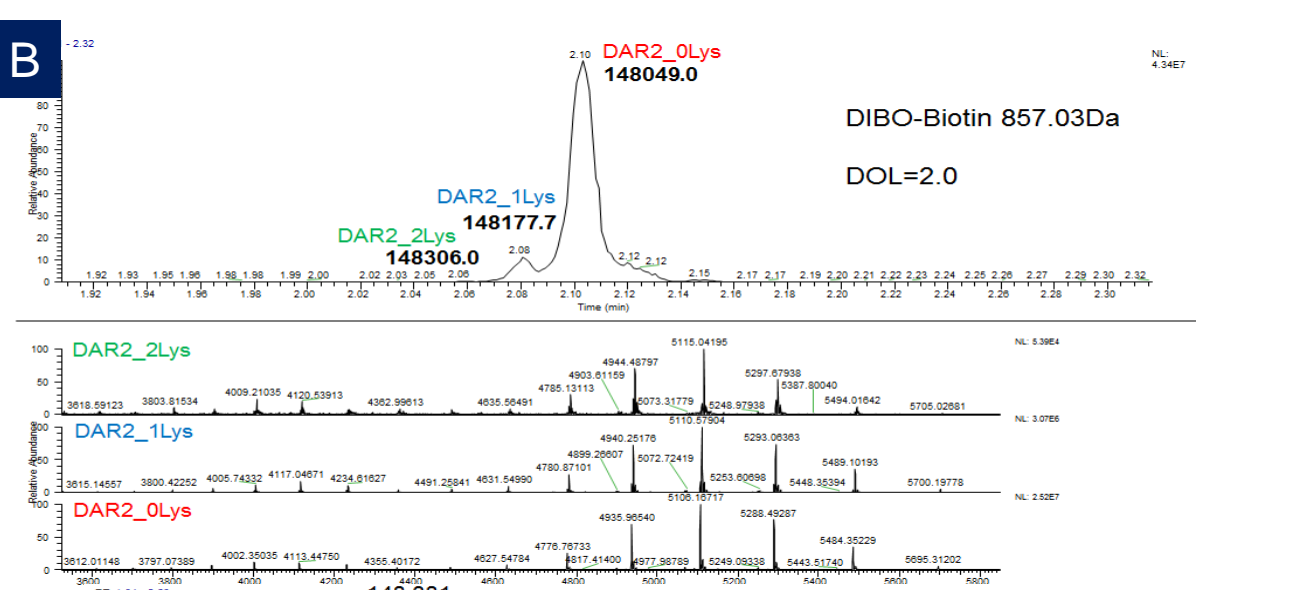
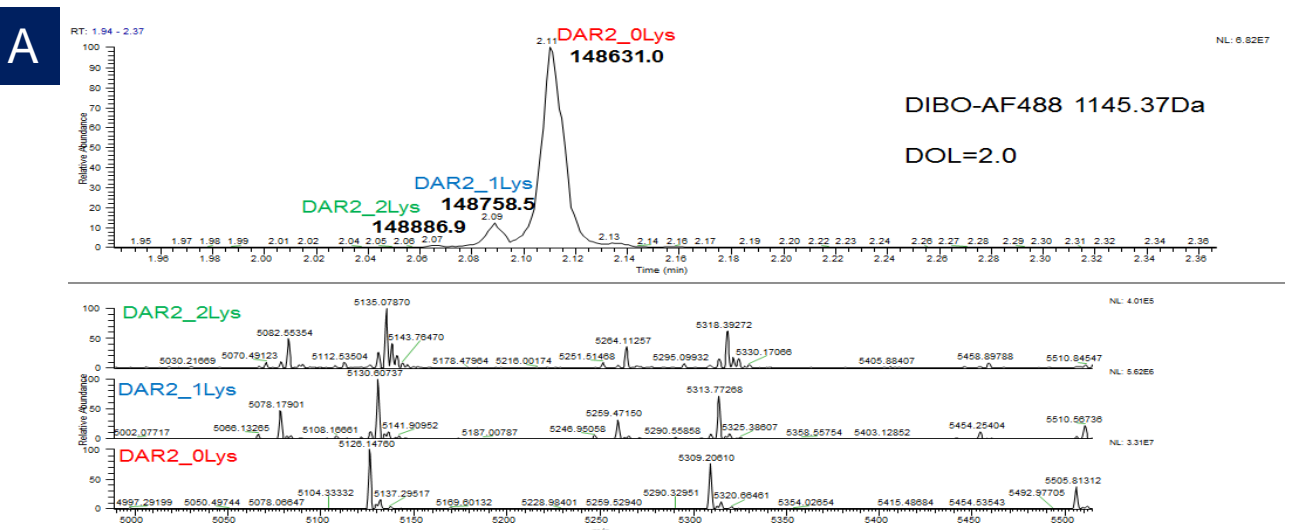


Figure 8. CEMS of NIST derived ADCs with A) AFF88, B) Biotin and C) dylight800 as payload on both N-glycan chains. CEMS separation conditions identical to conditions used for NIST mAb.

## CONCLUSIONS

- NIST mAb and NIST mAb derived ADC samples were separated and characterized using microfabricated glass chips in a CEMS experiment in less than 3 minutes by coupling the 908 ZipChip with Thermo Q Exactive HF MS with Biopharma option. No sample preparation and desalting was necessary.
- Charge variants based on different lysine content of the NIST mAb can be separated on the chip based CE device. The additional positive charges added by lysine to the C-terminal of the heavy chain increases the mobility and leads to a clear distinction of the charge variants.
- Most of the proteoforms listed in the NIST mAb Reference Material can be identified in this study, among them 5 glycosylated variants including G0F/G0F, G0F/G1F, G1F/G1F, G1F/G2F, and G2F/G2F from 2-lys, 1-lys, and 0-lys migration time frames.
- The relative sample amounts of the detected charge variants covers 3 orders of dynamic range.
- Despite the acidic separation condition, the NIST mAb exhibits a charge state envelope of 26-31, very close to native conditions.
- The NIST mAb derived ADCs exhibit as expected the same charge heterogeneity.
- The observed MW for the AF488 and biotin ADCs confirms one added payload per N-linked glycan.
- For the AF488 and Biotin as payloads, only the variant with 2 payloads were observed confirming the high yield of the click chemistry.
- The dylight800 variant shows a more complex pattern making further analysis necessary.

## REFERENCES

- John E. Schiel, Darryl L. Davis, and Oleg V. Borisov, "State-of-the-Art and Emerging technologies for Therapeutic Monoclonal Antibody Characterization: The NISTmAb Case Study", American Chemical Society (2015)
- Erin A. Redman, J. Scott Mellors, Jason A. Starkey, and Michael Ramsey, "Characterization of Intact Antibody Drug Conjugate Variants Using Microfluidic Capillary Electrophoresis – Mass Spectrometry" Anal. Chem. 88, 2220-2226 (2016)

## ACKNOWLEDGEMENTS

The authors gratefully acknowledge the help and support of Jennifer N. Sutton and Michael A. Blank from Thermo Fisher Scientific.

## TRADEMARKS/LICENSING

© 2017 Thermo Fisher Scientific Inc. All rights reserved. ZipChip is a trademark of 908 Devices 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.