

Evaluation of a microfluidic electrophoresis device coupled to an Orbitrap mass spectrometer for the characterization of biotherapeutic proteins.

Stephane Houel¹, Terry Zhang¹, Erin Redman², Scott Mellors², Aaron Bailey¹, Chris Petty², Aran Paulus¹, Kai Zhou¹, Jonathan Josephs¹,
¹Thermo Fisher Scientific, San Jose, CA, USA, 95134; ²908 Devices, Boston, MA, USA

ABSTRACT

Purpose: To evaluate the performance of a chip-based electrophoresis device coupled to an Orbitrap mass spectrometer for intact and sub-unit mass analyses and peptide mapping.

Results: Using the ZipChip device, the lysine variants of the NIST mAb were successfully identified at the intact and sub-unit level. For the glycoforms G0F/G1F and G1F/G1F of trastuzumab emtansine, the average DAR values were respectively 3.46 and 3.47, which are in accordance with previously published data. Using the ZipChip device, a sequence coverage in excess of 97% was observed for the light and heavy chains after a 10 min peptide mapping experiment.

INTRODUCTION

The discovery and development of biotherapeutics continues to accelerate. The complexity of these agents and the increasing requirements to characterize them for both safety and efficacy places a large burden on the analytical scientists tasked with these challenging demands. Intact and sub-unit mass analysis as well as peptide mapping are widely used to get insights on biotherapeutics. Often these assays are LC-MS based, but the development of a new microfluidic capillary electrophoresis device could offer a fast and sensitive orthogonal mode of separation. Here we evaluate the performance of a chip-based electrophoresis device coupled to an Orbitrap mass spectrometer for some of the major workflows used to characterize biotherapeutics

MATERIALS AND METHODS

Sample Preparation

The analyzed samples were the NIST mAb and trastuzumab emtansine. For intact mass analysis, the sample was simply diluted in water prior to analysis.

For sub-unit mass analysis, the sample was first diluted in Tris HCl 0.1M and digested with the IdeS protease. Subsequently, the sample was denatured with Guanidine 8M and reduced with DTT. The final step is a buffer exchange with the dilution buffer provided by 908 Devices in the peptide kit.

For peptide mapping, NIST mAb was denatured in guanidine HCl and Tris followed by reduction and alkylation with DTT/IAA. The alkylation was quenched with DDT following buffer exchange in 50 mM Tris using BioSpin™ 6 columns (Bio-Rad Laboratories). Trypsinization was performed at 37 °C for 30 min and quenched by lowering the pH with formic acid. Finally, the sample was diluted with the peptide dilution buffer provided in the 908 devices peptide kit.

Methods

Microfluidic electrophoresis

The ZipChip™ HR chip from 908 Devices (Boston, MA) was used for all experiments. For intact and subunit mass analysis and peptide mapping, field strengths of 500V, 220V and 400V were respectively used.

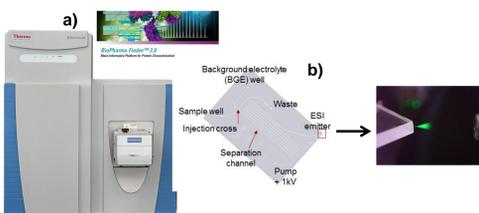
Mass Spectrometry

Mass spectrometer: A Thermo Scientific™ Q Exactive™ HF MS operated in high mass range (HMR) mode was used for the intact mass analysis and in standard mode operation for peptide mapping. A Thermo Scientific™ Q Exactive™ Plus in protein mode was used for the sub-unit mass analysis.

Data Analysis

For all of the experiments, Thermo Scientific™ BioPharma Finder™ 2.0 software was used.

Figure 1. a) Data were collected using the ZipChip device coupled to a Q Exactive Plus or Q Exactive HF instrument and processed with BioPharma Finder software. **b)** The ZipChip HR is a twenty two centimeter etched channel with a neutral hydrophilic coating with an integrated nano-electrospray emitter



RESULTS

A) Intact Mass Analysis

1) NIST mAb

Figure 2. a) Electropherogram of NIST mAb (0.1 ug/uL). **b)** Spectra of the NIST mAb and the lysine variants. **c)** Zoom in the region of the charge +30 ions.

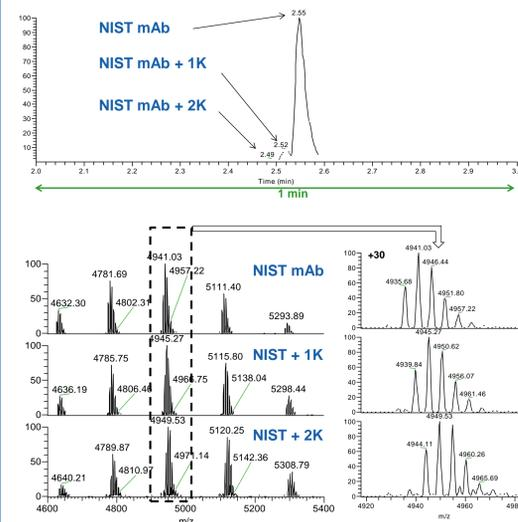


Table 1. Deconvoluted masses of the NIST mAb and lysine variants major glycoforms.

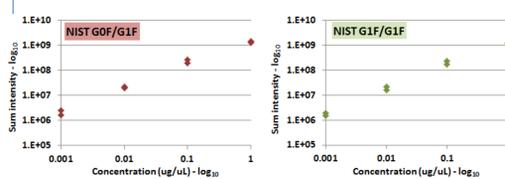
Protein Name	Modification	Average Mass	Theoretical Mass (Da)	Matched Mass Error (ppm)	Sum Intensity	Relative Abundance	Fractional Abundance
NIST		148038.8	148037.1	11.5	1.28E+09	58.82	16.61
NIST_plus1K	G0F/G0F	148155.6	148155.3	1.9	7.66E+07	3.51	0.99
NIST_plus2K		148295.4	148293.5	13.1	1.04E+07	0.48	0.14
NIST		148200.0	148199.3	5.0	2.18E+09	100.00	28.24
NIST_plus1K	G0F/G1F	148327.4	148327.4	0.1	1.39E+08	6.36	1.80
NIST_plus2K		148456.9	148455.6	9.0	1.56E+07	0.72	0.20
NIST		148362.1	148361.2	6.0	1.82E+09	83.26	23.51
NIST_plus1K	G1F/G1F	148489.7	148489.4	2.5	1.18E+08	5.40	1.53
NIST_plus2K		148618.4	148617.5	5.6	1.51E+07	0.69	0.20
NIST		148522.7	148523.5	5.8	1.00E+09	46.00	12.99
NIST_plus1K	G1F/G2F	148652.3	148651.7	3.6	5.97E+07	2.74	0.77
NIST_plus2K		148779.4	148779.9	3.1	8.07E+06	0.37	0.10
NIST		148684.6	148685.7	7.2	4.89E+08	22.21	6.27
NIST_plus1K	G2F/G2F	148811.9	148813.9	13.4	2.90E+07	1.33	0.38
NIST_plus2K		148942.0	148942.0	0.4	3.49E+06	0.16	0.05

0.65 nL of the NIST mAb at 0.1 ug/uL was injected and the three lysine variants are almost baseline separated. The shift in mass can be observed in the zoom in the region of the +30 charge state ions. The fast and almost base line separation of the lysine variants using the ZipChip combined with the high mass accuracy of the Orbitrap analyzer translated to the identification of 15 different glycoforms with a mass error below 14 ppm.

Table 2. Deconvoluted masses and intensities of the NIST mAb for the glycoforms G0F/G1F and G1F/G1F at different concentrations.

Concentration (ug/uL)	Injection Volume = 0.65 nL				Injection Volume = 0.65 nL				
	Average Mass (Da)	Theoretical Mass (Da)	Matched Mass Error (ppm)	Sum Intensity	Average Mass (Da)	Theoretical Mass (Da)	Matched Mass Error (ppm)	Sum Intensity	
0.001	148199.55	148199.26	2.0	1.66E+06	0.001	148362.39	148361.20	8.0	1.55E+06
0.01	148199.59	148199.26	2.2	2.26E+07	0.01	148362.02	148361.20	5.5	2.28E+07
0.1	148199.47	148199.26	1.5	1.95E+07	0.1	148361.14	148361.20	0.4	1.63E+07
1	148200.60	148199.26	9.1	2.70E+08	1	148361.93	148361.20	4.9	2.35E+08
	148200.36	148199.26	7.4	1.96E+08		148362.41	148361.20	8.2	1.71E+08
1	148199.01	148199.26	1.7	1.43E+09	1	148362.74	148361.20	10.4	1.18E+09
	148199.55	148199.26	4.1	1.29E+09		148360.64	148361.20	3.8	1.20E+09

Figure 3. graphs representing the sum of the intensity of the deconvoluted masses versus the concentration for the NIST mAb glycoforms G0F/G1F and G1F/G1F.



The low sample consumption and high sensitivity of the intact mass analysis assay with the ZipChip are illustrated with the dilution series of the NIST mAb. Even at 0.001 ug/uL when only 0.65 nL are injected, glycoforms G0F/G1F and G1F/G1F are identified at less than 10 ppm.

2) Trastuzumab Emtansine

Figure 3. Data visualization in BioPharma Finder 2.0 software of the lysine-linked ADC.

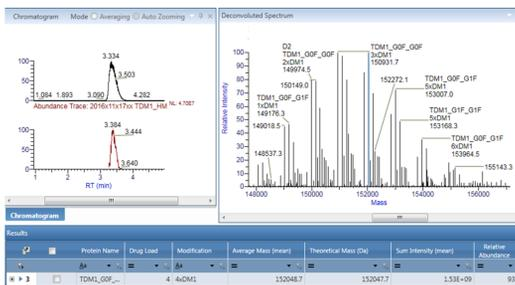


Table 3. Average DAR calculation for glycoforms G0F/G1F and G1F/G1F of trastuzumab emtansine

Sample Name	Sample Concentration (ug/uL)	Average Mass (Da)	Theoretical Mass (Da)	Matched Mass Error (ppm)	Sum Intensity	Relative Abundance	Fractional Abundance
Trastuzumab Emtansine	1.0	148794.5	148794.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	148918.5	148918.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	149042.5	149042.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	149166.5	149166.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	149290.5	149290.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	149414.5	149414.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	149538.5	149538.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	149662.5	149662.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	149786.5	149786.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	149910.5	149910.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	150034.5	150034.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	150158.5	150158.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	150282.5	150282.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	150406.5	150406.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	150530.5	150530.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	150654.5	150654.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	150778.5	150778.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	150902.5	150902.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	151026.5	151026.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	151150.5	151150.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	151274.5	151274.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	151398.5	151398.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	151522.5	151522.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	151646.5	151646.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	151770.5	151770.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	151894.5	151894.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	152018.5	152018.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	152142.5	152142.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	152266.5	152266.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	152390.5	152390.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	152514.5	152514.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	152638.5	152638.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	152762.5	152762.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	152886.5	152886.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	153010.5	153010.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	153134.5	153134.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	153258.5	153258.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	153382.5	153382.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	153506.5	153506.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	153630.5	153630.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	153754.5	153754.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	153878.5	153878.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	154002.5	154002.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	154126.5	154126.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	154250.5	154250.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	154374.5	154374.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	154498.5	154498.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	154622.5	154622.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	154746.5	154746.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	154870.5	154870.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	154994.5	154994.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	155118.5	155118.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	155242.5	155242.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	155366.5	155366.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	155490.5	155490.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	155614.5	155614.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	155738.5	155738.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	155862.5	155862.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	155986.5	155986.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	156110.5	156110.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	156234.5	156234.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	156358.5	156358.5	0.0	1.00E+09	100.00	100.00
Trastuzumab Emtansine	1.0	156482.5	156				