

Poster Reprint

**ASMS 2024**  
**Poster number MP 665**

# Peptide Mapping of Tryptic Digests for mAbs using a novel ECD cell on the 6545XT AdvanceBio LC/Q-TOF Mass Spectrometer

Stephen Sciuto<sup>1</sup>, Maozi Liu<sup>1</sup>, Rachel Franklin<sup>2</sup>, Jerry Han<sup>1</sup>

<sup>1</sup>Agilent Technologies, Inc., Santa Clara, CA

<sup>2</sup>Agilent Technologies, Corvallis, Oregon

## Introduction

Post-translation modifications (PTMs) on monoclonal antibodies (mAb) play an important role in the safety, efficacy, and binding of a therapeutic to its target [1]. Common examples of PTMs that seek to be identified include glycosylation and phosphorylation. One challenge in the identification of these PTMs is that dissociation techniques such as collision-induced dissociation (CID) can break apart these fragile modifications. Here, we describe the use of an electron-based dissociation technique (ExD) that supplies low-energy electrons for the fragmentation of tryptic digests that contain glycosylation at asparagine consensus sites (NXS/T). The fragmentation of trastuzumab tryptic digest on a 6545XT AdvanceBio LC/Q-TOF using CID is compared to illustrate that CID alone will fragment these labile PTMs.

## Experimental

Tryptic digests for mAbs were generated using the Agilent AssayMAP Bravo system using either an in-solution digestion or single-pot (SP3) protocol. Following digestion, peptides were reconstituted in 0.1% formic acid in water at a concentration of  $\sim 0.5$   $\mu\text{g}/\mu\text{L}$ . Peptides were added to a G7167B multisampler at 6°C inside a 1290 Infinity II LC system. LC/MS data was collected on a 6545XT equipped with an ExD cell and searched using a pre-released version of MassHunter BioConfirm 12.1 software. For tuning, the Extended Dynamic Range (2 GHz) mode was used with the 100 – 3000  $m/z$  range.

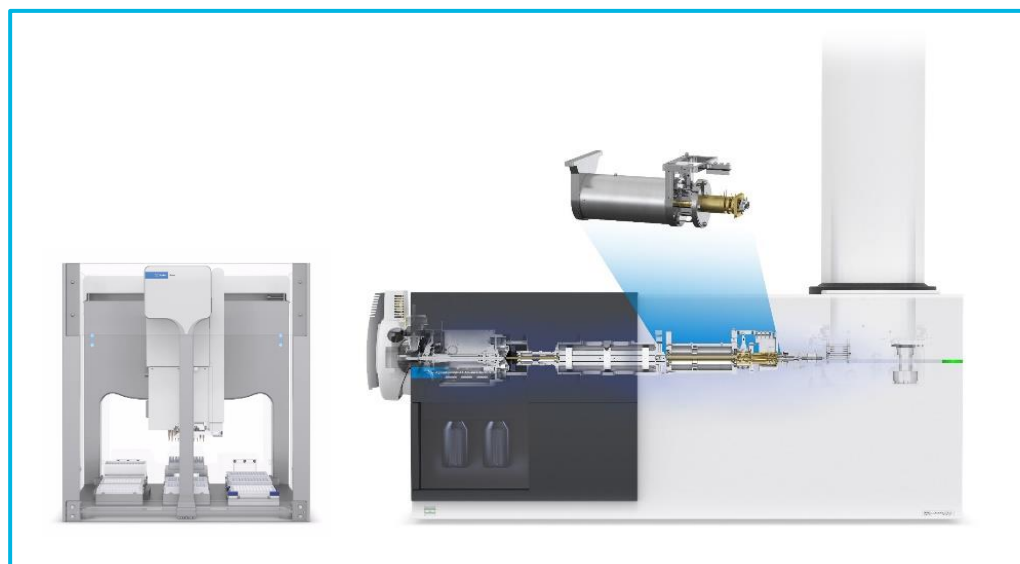


Figure 1. AssayMAP Bravo (left) and Agilent 6545XT AdvanceBio LC/Q-TOF with ExD cell (right).

## Experimental

### LC Conditions

Solvent A	Water with 0.1% FA
Solvent B	ACN with 0.1% FA
Gradient	0-90 min, 2-40% B 90.5-94 min, 80% B 94.5-100 min, 2-40% B 100-104 min, 80% B
Injection volume	6 $\mu\text{L}$
Flow rate	0.3 mL/min
Column temperature	40°C

### MS Conditions

Gas temperature	250 °C
Drying gas	10 L/min
Nebulizer	25 psi
Sheath gas temperature	250 °C
Sheath gas flow	12 L/min
Vcap	3500 V
Nozzle voltage	0 V
Fragmentor	170 V
Skimmer	65 V
Reference mass	922.0098
MS <sup>1</sup> range	200 – 3000 $m/z$
Acquisition rate	3 spectra/sec
MS/MS range	100 – 3000 $m/z$
Acquisition rate	3 spectra/sec
Isolation width	Medium ( $\sim 4$ $m/z$ )
Precursors/cycle	Top 5
Threshold for MS/MS	3000 counts and 0.001%
Precursor charge	2+, 3+, >3+
Target	50,000 counts/spectrum

### Glycopeptide signal in base peak chromatogram

Glycopeptides for NIST and trastuzumab typically contribute <10% relative contribution to the base peak chromatogram on the Agilent 6545XT AdvanceBio LC/Q-TOF with ExD cell. However, these signals still allow for excellent sequence coverage along the peptide backbone for these labile molecules (see Figure 3 and 4).

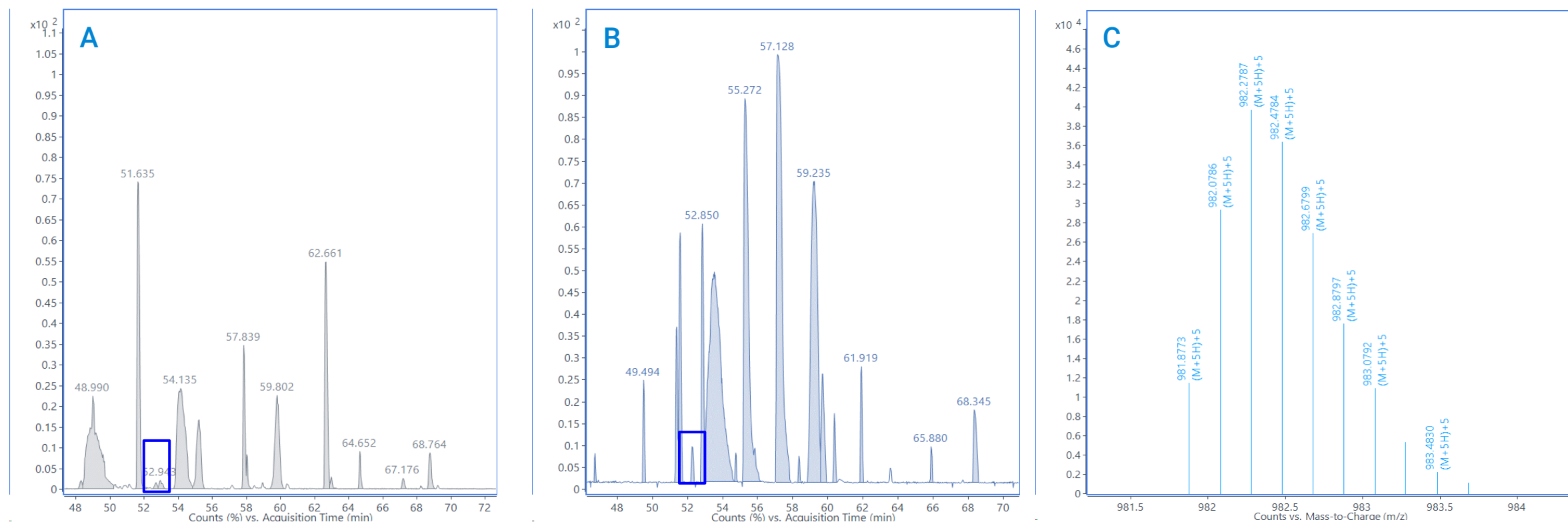


Figure 2. Zoomed-in region of the base peak chromatogram illustrating the relative abundance of the glycopeptide signal (blue box) for (A) NIST and (B) Trastuzumab. A representative MS<sup>1</sup> spectrum is displayed for Trastuzumab in (C) in centroid mode. MS/MS spectra for these glycopeptides are measured using ExD or CID mode with the MS criteria displayed in Table 1.

### NIST glycopeptide MS/MS on 6545XT with ExD cell

Tryptic digests of NIST mAb in ExD mode resulted in a sequence coverage of ~99%. In addition, glycopeptide MS/MS illustrated excellent sequence coverage along the peptide backbone (Figure 3). Noteworthy is the appearance of c-type peptide fragments that contain G1F (solid blue box) and those without G1F (dashed blue box).

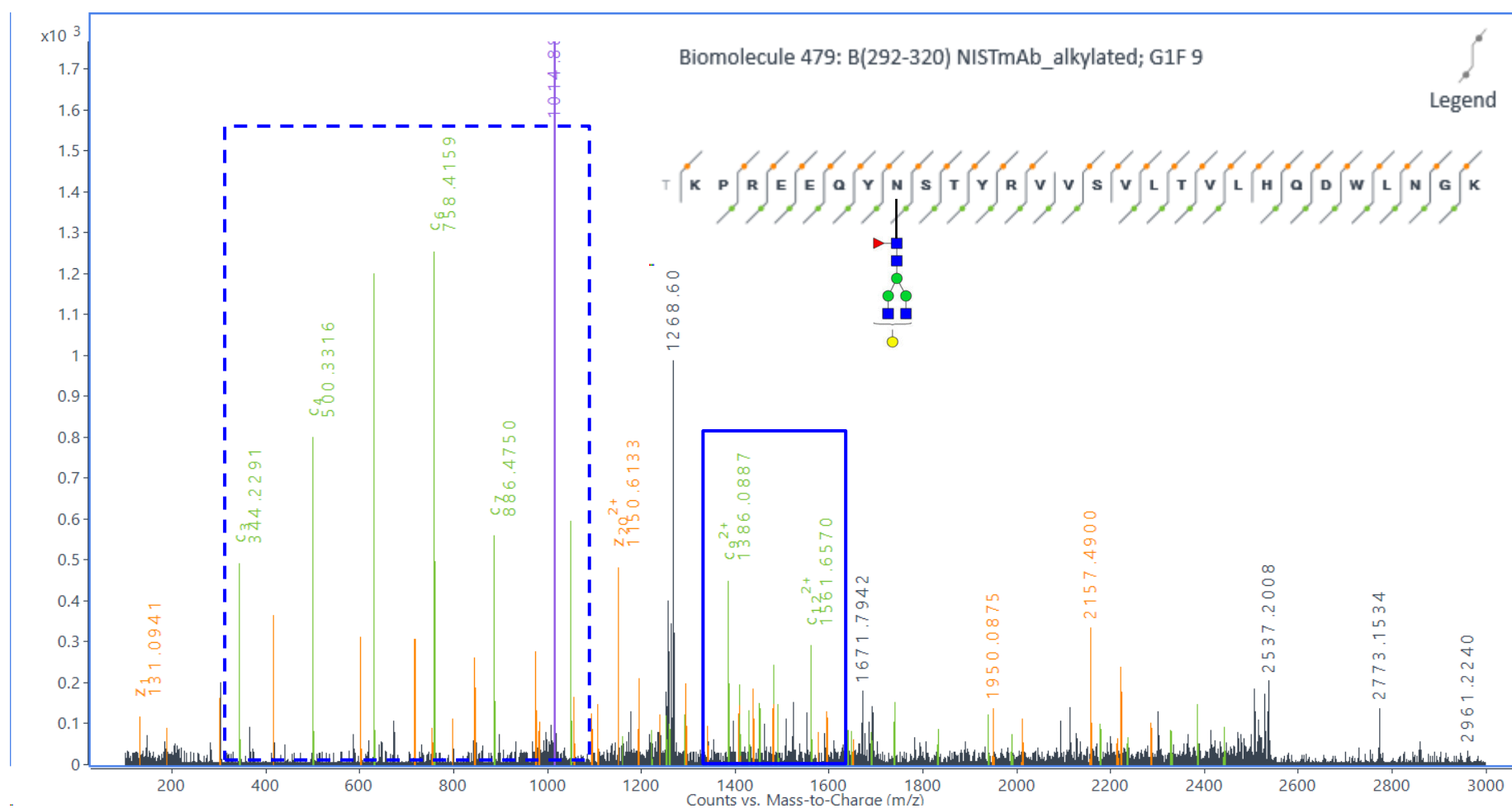


Figure 3. MS/MS spectrum and peptide ladder for TKPREEQYNSTYRVVSVLTVLHQDWLNGK with G1F modification.

## Trastuzumab glycopeptide MS/MS on 6545XT with ExD cell

Tryptic digests of trastuzumab in ExD mode resulted in a sequence coverage of ~99%. Glycopeptide MS/MS illustrated excellent sequence coverage along the peptide backbone (Figure 4A), though CID generated the HexNAc glycan fragment at  $m/z$  204.086 (Figure 4B). Like NIST, there is an abundance of c-type fragment ions that contain the glycopeptide. In addition, there are z-type fragment ions with (blue box) and without the glycopeptide (blue dashed box). A zoomed-in region also shows the mass difference between  $z_{21}^{2+}$  (with GOF) and  $z_{20}^{2+}$  (without GOF) fragment ions.

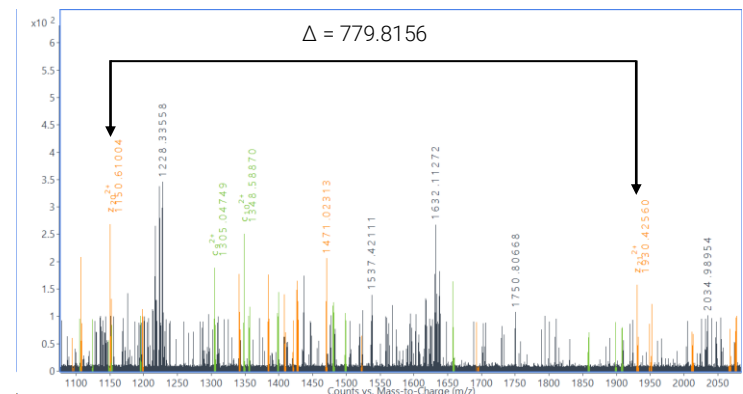
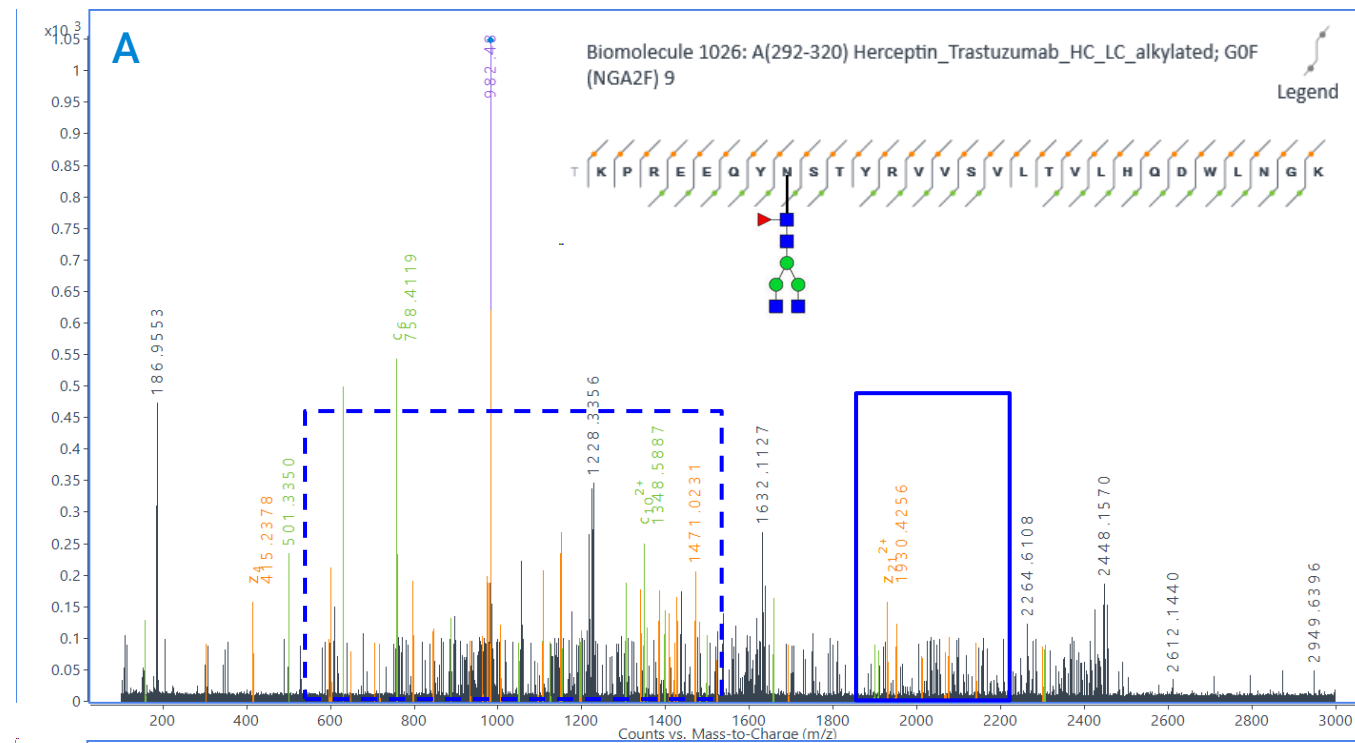


Figure 4. (A) MS/MS spectrum and peptide ladder for TKPREEQYNSTYRVVSVLTVLHQDWLNGK with GOF modification in the 5+ charge state operated in ExD mode. The solid blue box indicates fragment ions with GOF attached whereas the dashed blue box shows fragment ions without GOF. The zoomed-in region above shows  $z_{21}^{2+}$  (with GOF) and  $z_{20}^{2+}$  (without GOF).

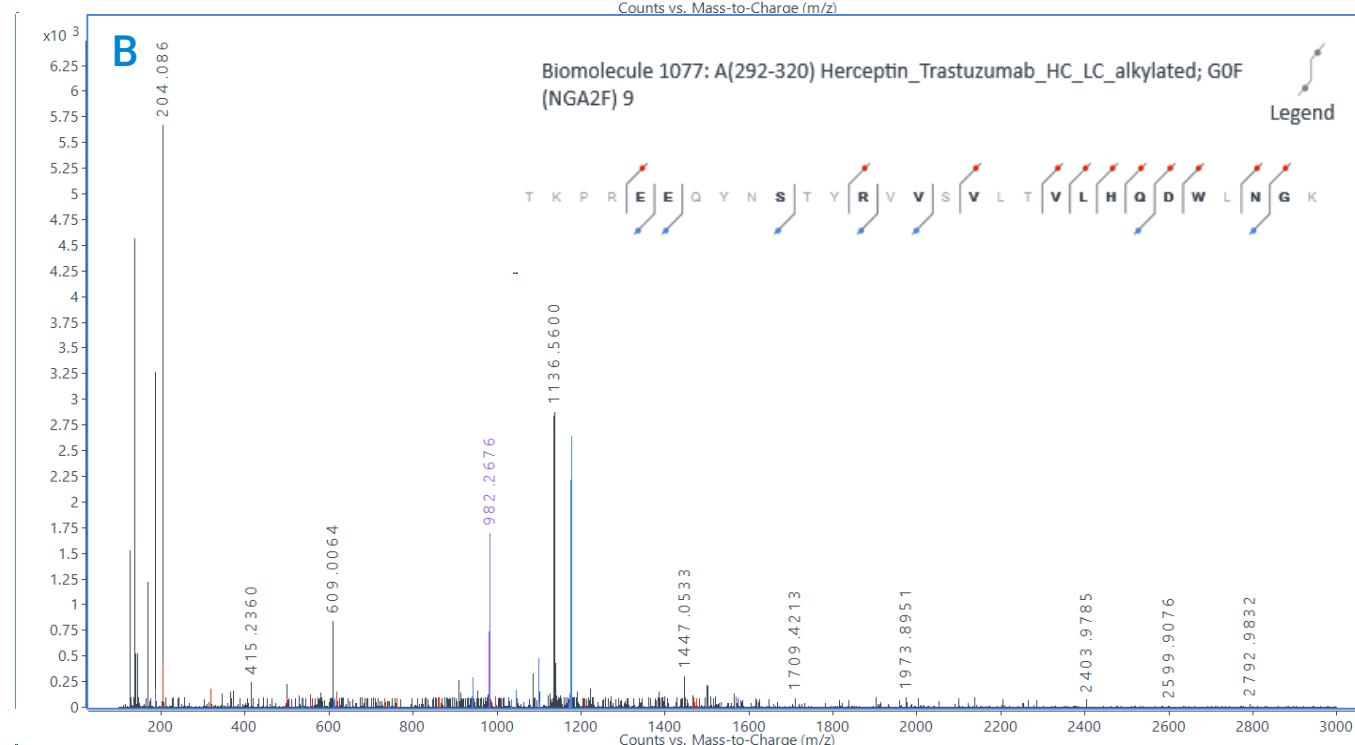


Figure 4. (B) MS/MS spectrum and peptide ladder for TKPREEQYNSTYRVVSVLTVLHQDWLNGK with GOF modification in the 5+ charge state operated in CID mode. Fragmentation of this glycopeptide generates the HexNAc glycan fragment at  $m/z$  204.086 with fewer fragments along the peptide backbone.

## Conclusions

### The Agilent 6545XT AdvanceBio LC/Q-TOF with ExD cell enables glycopeptide analysis with excellent sequence coverage assessed using MassHunter BioConfirm 12.1

- LC/MS conditions were optimized to detect low-abundant glycopeptide species from NIST and trastuzumab
- MS/MS spectra for glycopeptides using ExD show peptide fragments with intact glycans and excellent sequence coverage along the peptide backbone
- For CID mode, no intact glycan fragments are observed. Instead, HexNAc glycan fragment at  $m/z$  204.086 is observed

## References

- [1] Zheng K, Bantog C, Bayer R. The impact of glycosylation on monoclonal antibody conformation and stability. *MAbs*. 2011 Nov-Dec;3(6):568-76. doi: 10.4161/mabs.3.6.17922. Epub 2011 Nov 1. PMID: 22123061; PMCID: PMC3242843.

<https://www.agilent.com/en/promotions/asms>

This information is subject to change without notice.

DE29227673