

# Glycoprotein characterization combining native Mass Spectrometry and Direct Mass Technology

Weijing Liu, Ping Yip, Christopher Mullen, Claire Dauly, Rosa Viner  
Thermo Fisher Scientific, San Jose, CA, United States, 95134

## Abstract

**Purpose:** Using proton transfer charge reduction, hybrid fragmentation, and Direct Mass Technology to unravel glycoprotein complexity

- Methods:**
1. Proton transfer charge reduction (PTCR)
  2. Electron transfer dissociation supplemented with HCD
  3. Direct Mass Technology mode

**Results:** Complementary mass spectrometry analyses including Native, PTCR, Direct Mass Technology mode, and EThcD could comprehensively characterize heavily glycosylated proteins.

## Introduction

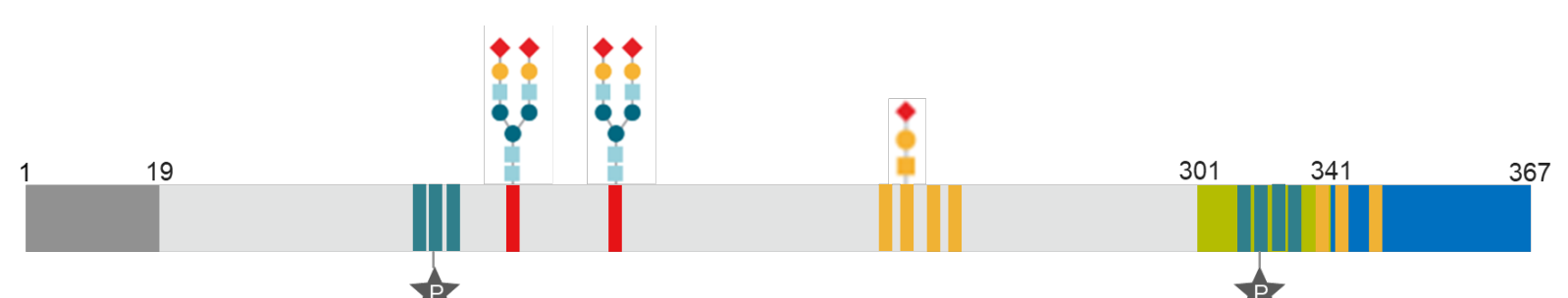
Glycoproteins are actively involved in physiological functions as well as certain diseases. However, glycoprotein characterization using native and native top-down mass spectrometry remains challenging as its heterogeneity leading to complex spectrum.

In this study, we evaluated a few novel techniques including Direct Mass Technology mode and proton transfer charge reduction (PTCR) to unravel proteoforms of Human Fetuin A, a heavily glycosylated and phosphorylated protein, comprehensively under native condition. We further expand the application to heavily glycosylated Spike protein variants for determining their MW and oligomeric states.

## Materials and methods

### Sample preparation

Human Fetuin A from human plasma was purchased from Sigma-Aldrich. Fetuin was buffer exchanged into ammonium acetate with Amicon 10K-MWCO or 30K-MWCO.



Average MW 37343.23 (w/o signal peptide and PTMs)

Signal Peptide: AA1-18

A-chain: AA19-300

Propeptide: AA301-340

B-chain: AA341-367

N-glycosylation: N156, and N176

Phosphorylation: S134, S135, S138, T319, S325, S328, S330

O-glycosylation: S256, T270, S280, S293, T339, T341, S346

Cys-Cys: 12 disulfide bonded cysteines

R340 can be missing

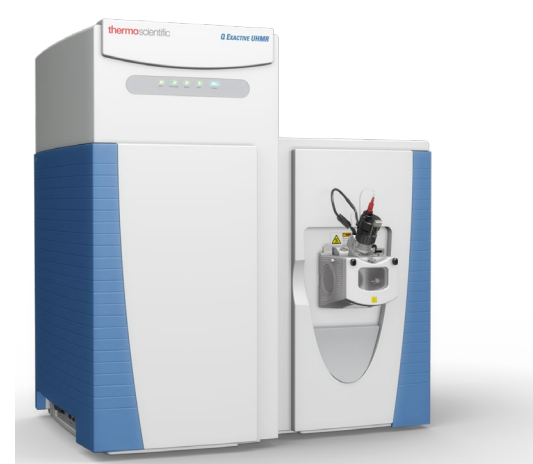
Spike protein samples expressed in HEK 293 cells were provided by Dr. Ganesh Anand, Penn State University.

### Test methods

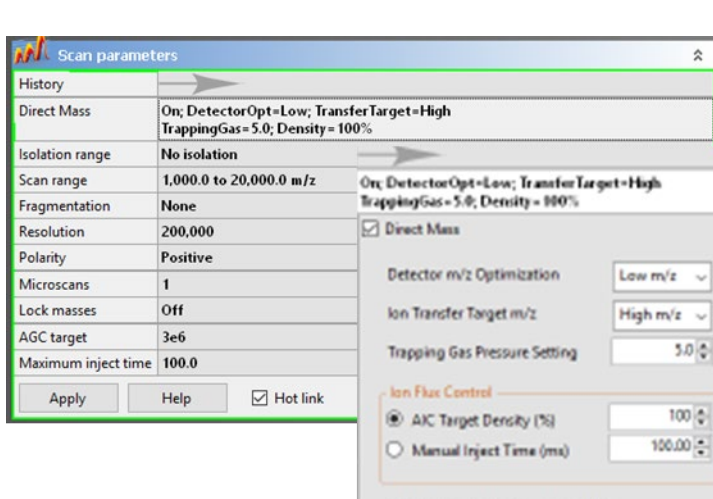
Native MS and Direct Mass Technology mode were performed on Thermo Scientific™ Q Exactive™ UHMR mass spectrometer. PTCR and EThcD analyses were performed on Thermo Scientific™ Orbitrap Eclipse™ mass spectrometer.



- Native MS
- Native top-down
- PTCR
- ETD
- UVPD

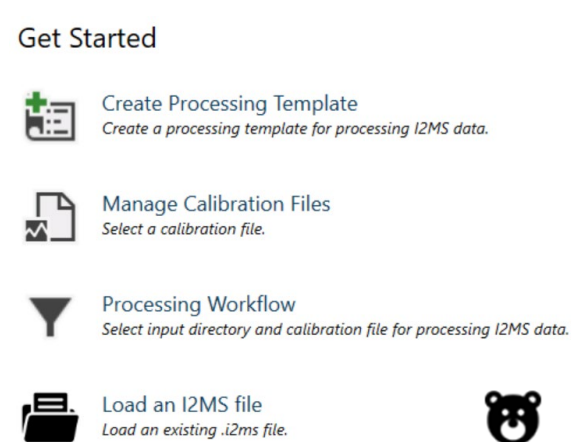


- Native MS
- Native top-down
- Direct Mass Technology



### Data Analysis

Data were analyzed using Thermo Scientific™ BioPharma Finder™ 4.1 and STORboard (Proteinaceous) software.



## Results

### Proton transfer charge reduction

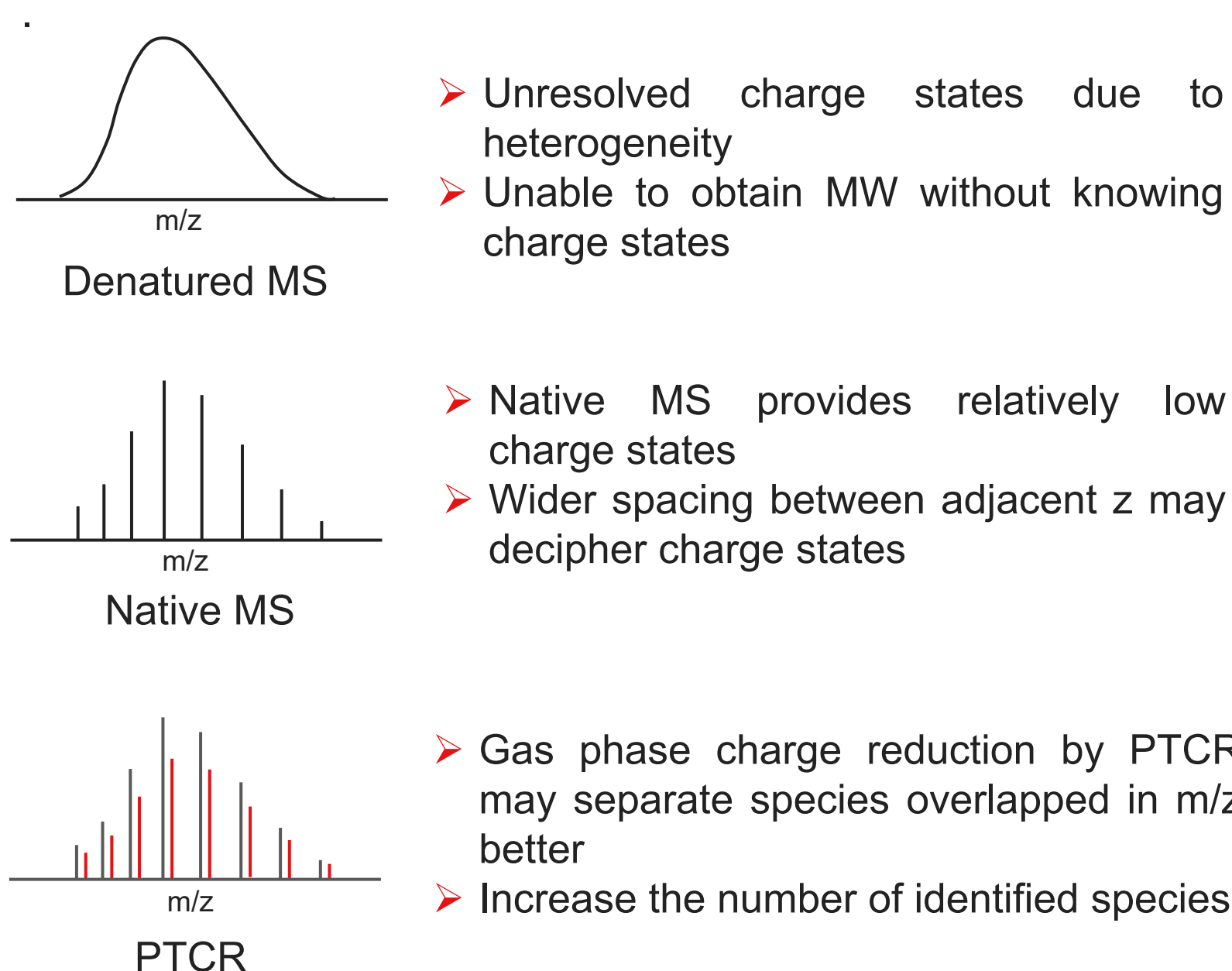
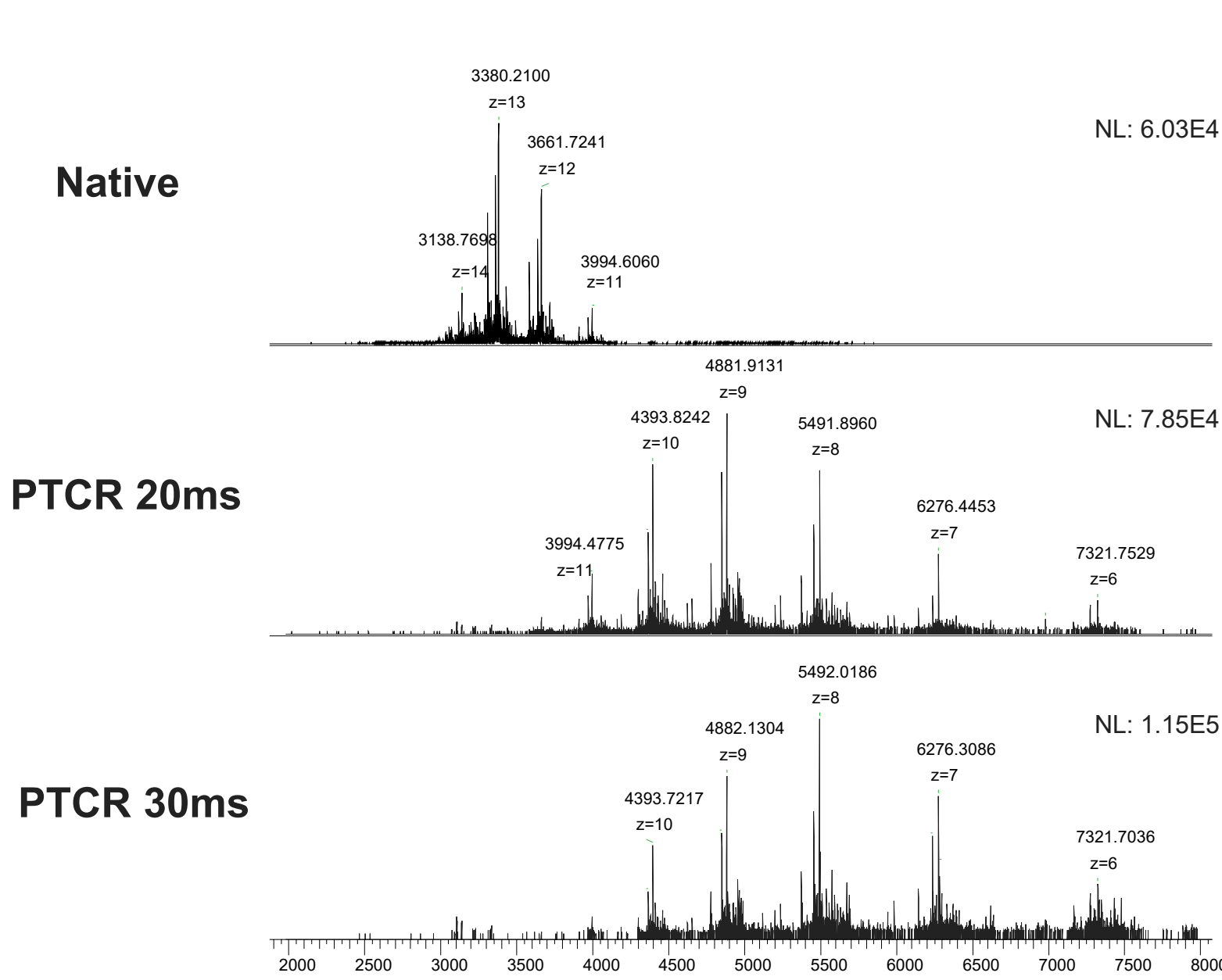
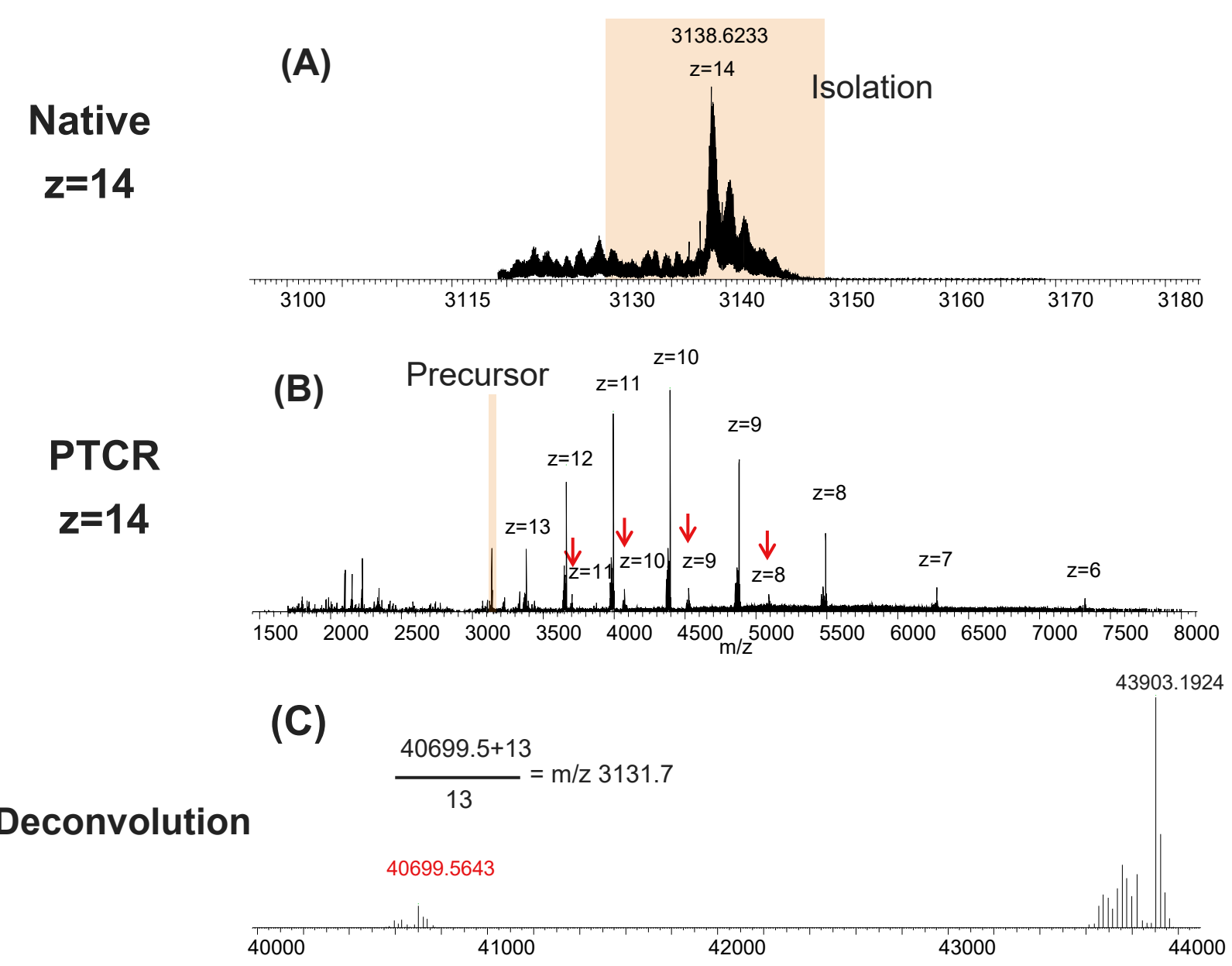


Figure 1. Comparison of native MS spectrum and PTCR MS spectra at 20 and 30 ms reaction time.



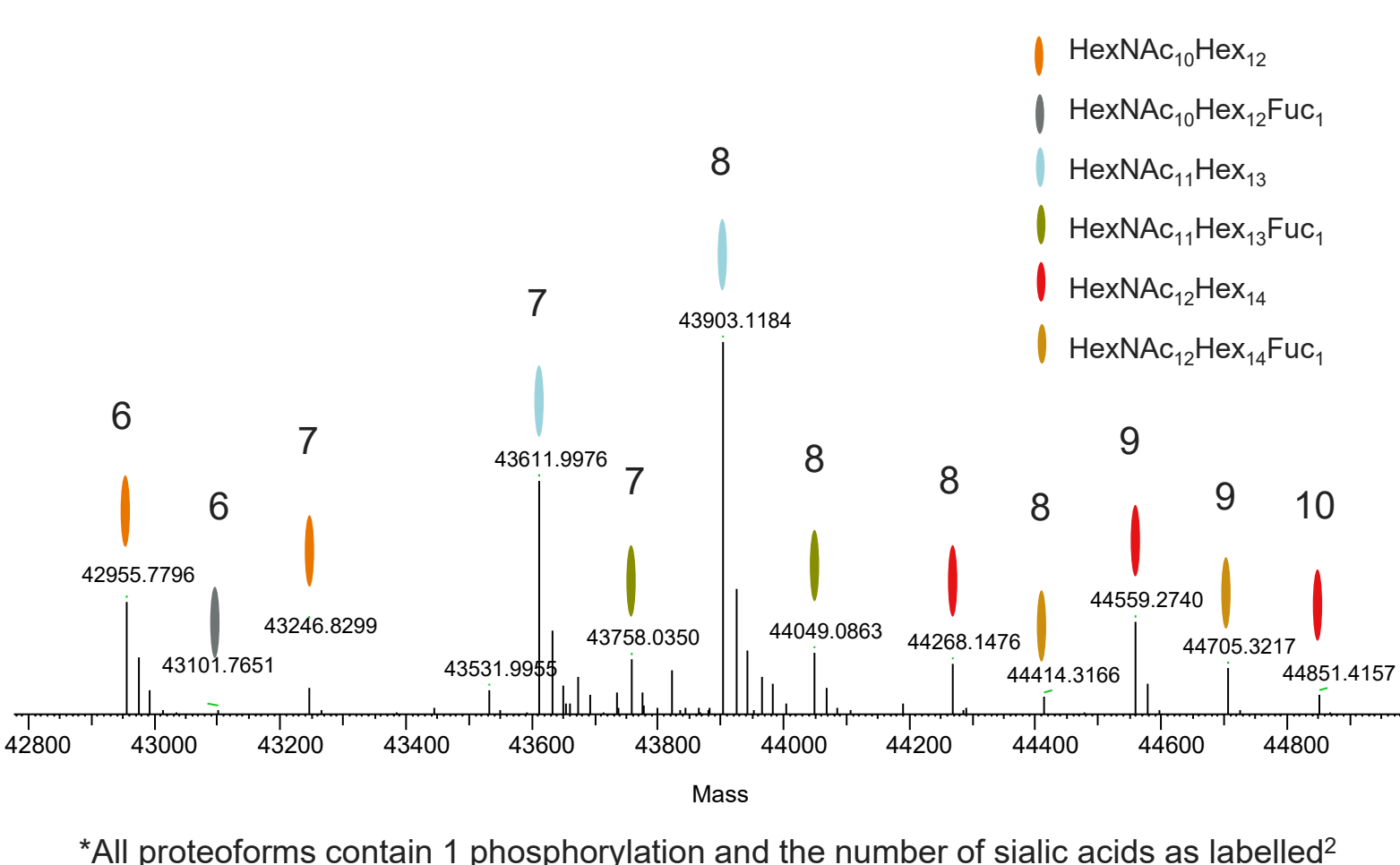
- PTCR<sup>1</sup> pushes the charge envelope to higher m/z and lower charge states via charge reduction
- Spectra reflects the sample complexity

Figure 2. (A) Zoom-in of z=14+ in native MS spectrum; (B) PTCR MS spectrum with 20 Th isolation width on z=14+; (C) Deconvoluted spectrum from PTCR scan.



- PTCR separates proteoforms previously overlapped in the full scan and thus increases number of identified glycoforms.

Figure 3. Fetuin A proteoforms identified using PTCR



- Accurate monoisotopic mass unambiguously deciphers different numbers of phosphorylation, HexNAc, Hex, and Fucose on each glycoform.

## Direct Mass Technology

- **Conventional MS:** Charge must be deduced from spacing between adjacent peaks of charge envelope or isotopic peaks
- **Challenges:** Heterogeneity causes loss of charge state resolution – particularly for large or heavily PTM molecules
- **Direct Mass Technology mode:** m/z and z are simultaneously measured for single ion

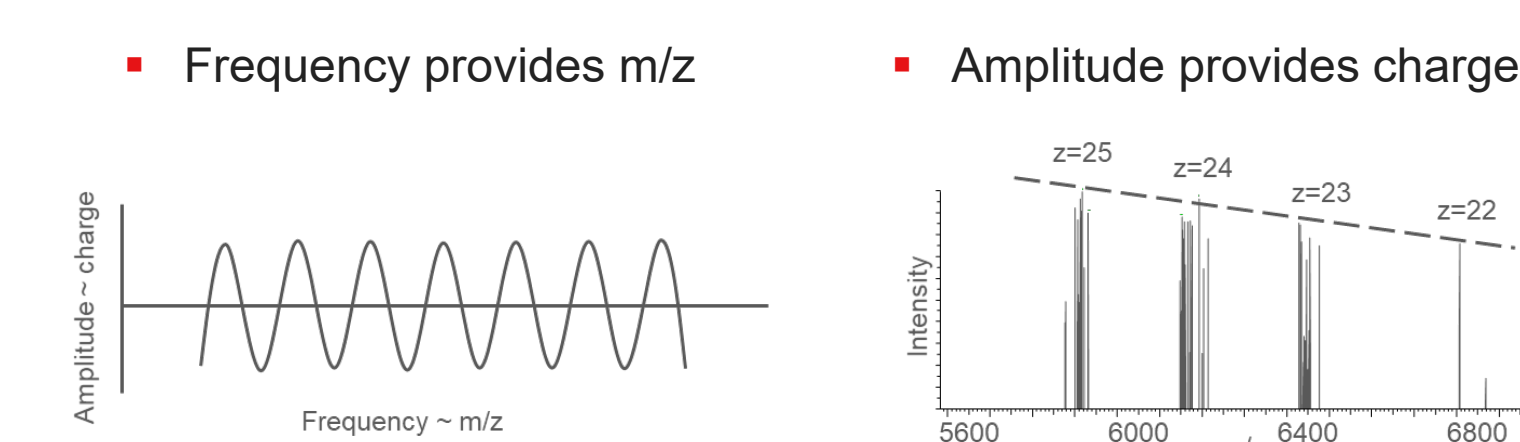
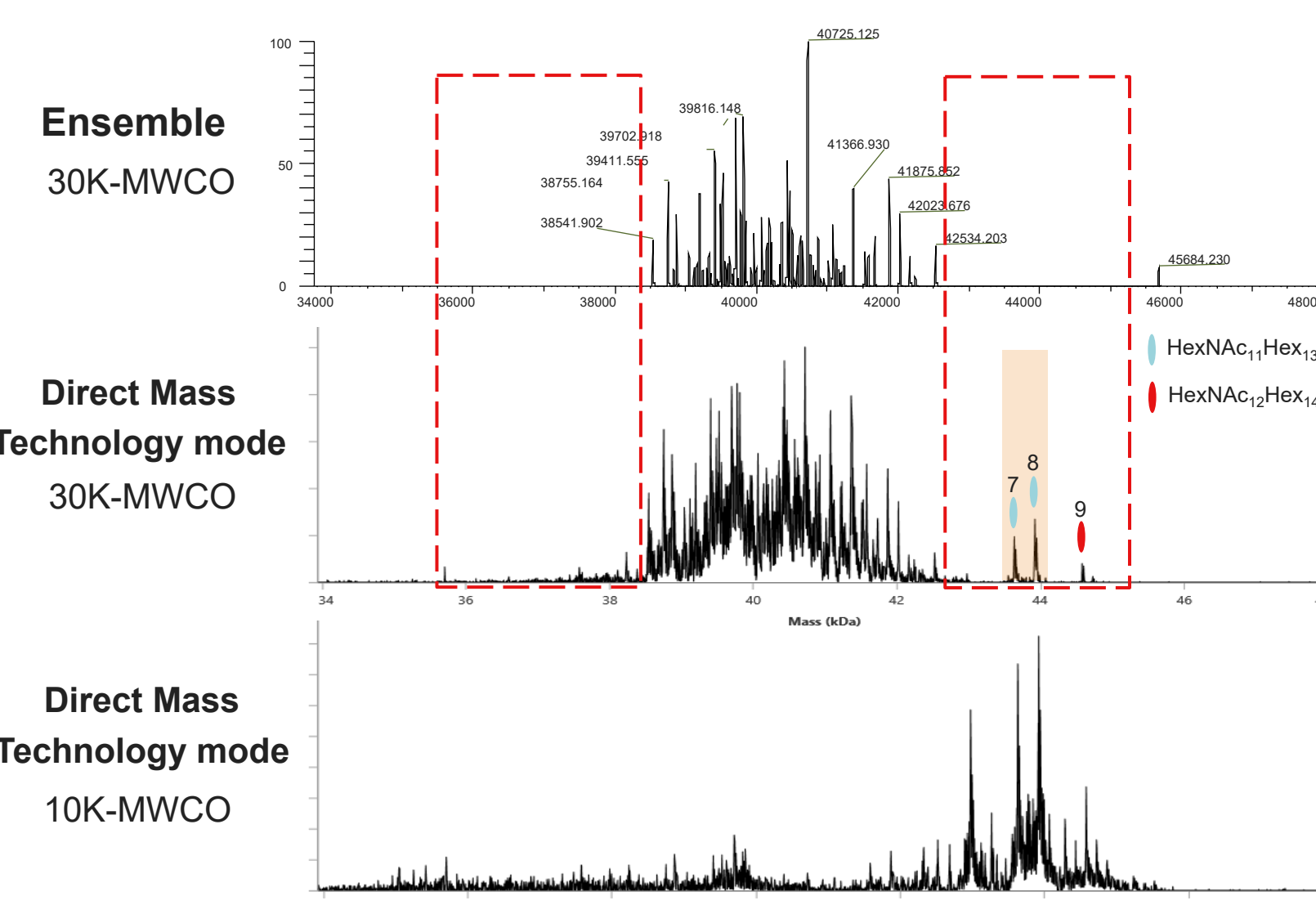
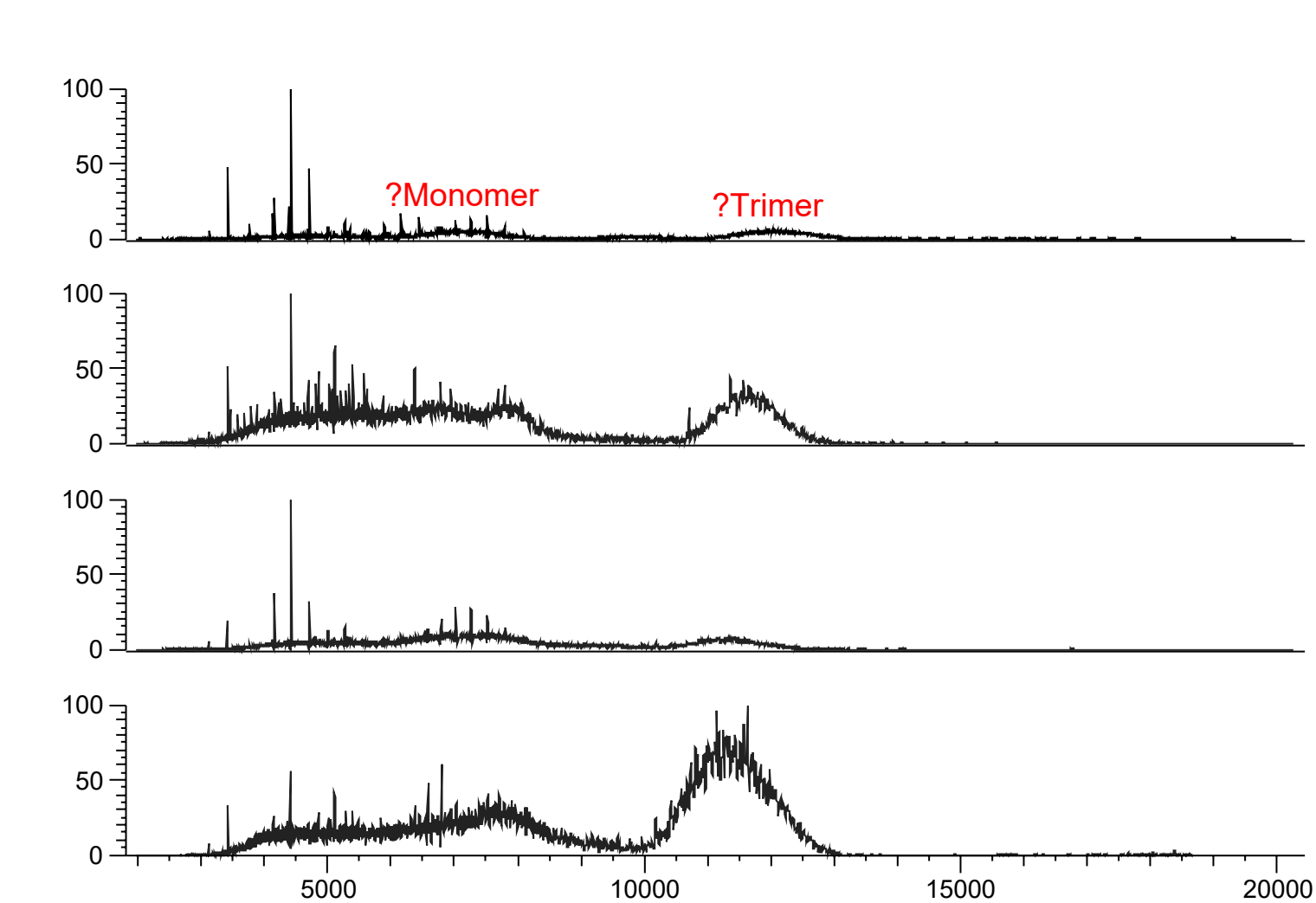


Figure 4. Comparison between spectra from ensemble measurement and Direct Mass Technology mode



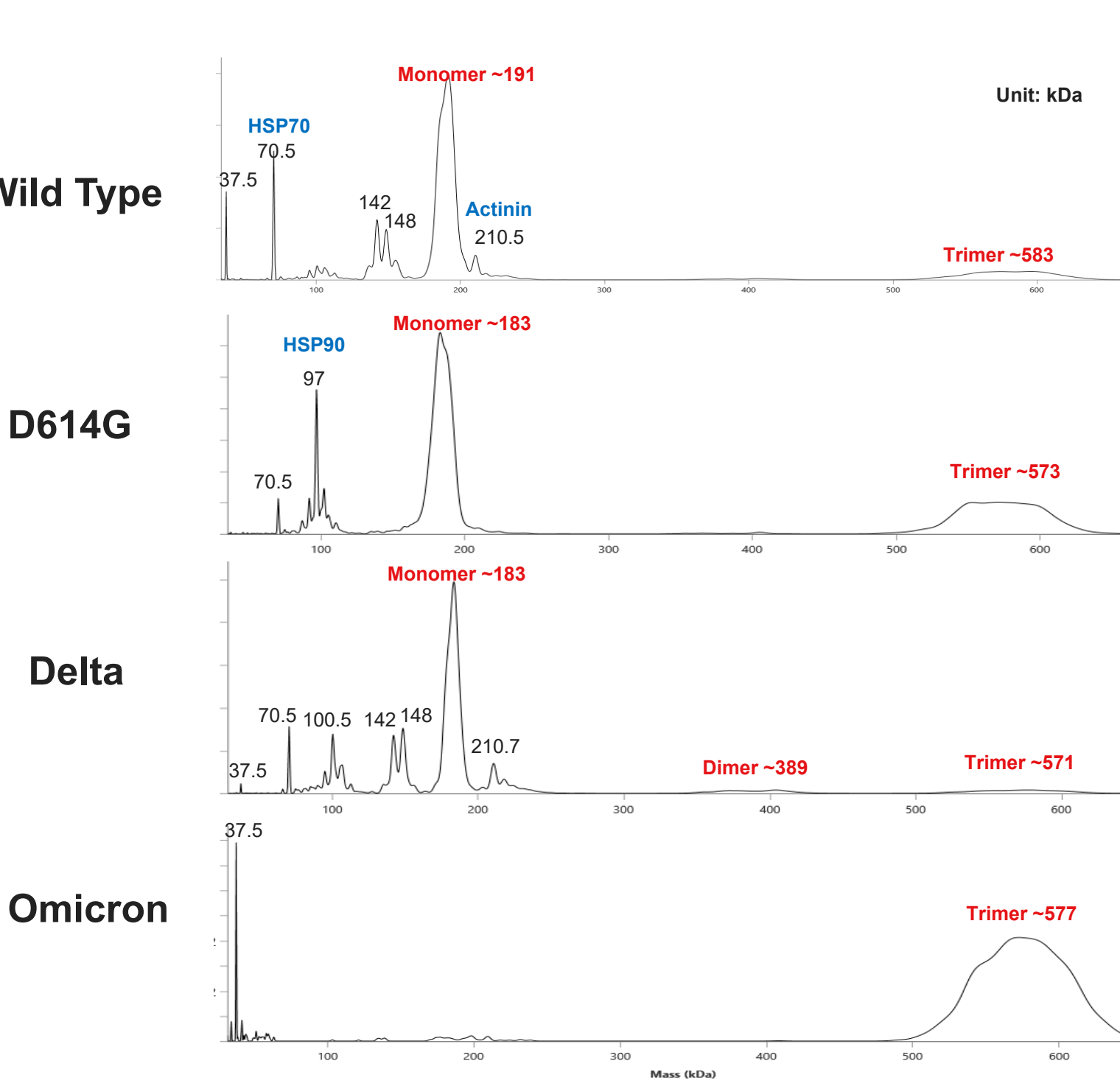
- Fetuin cleaned by 30K-MWCO indicate main glycoforms distributing from 38.5 kDa to 42.5 kDa
- Results demonstrate exceptional sensitivity and dynamic range of Direct Mass Technology mode

Figure 5. Ensemble MS Analysis of SARS-CoV-2 Spike Variants



- Ensemble measurement shows hump above and below m/z 10,000 potentially indicating trimer monomer respectively

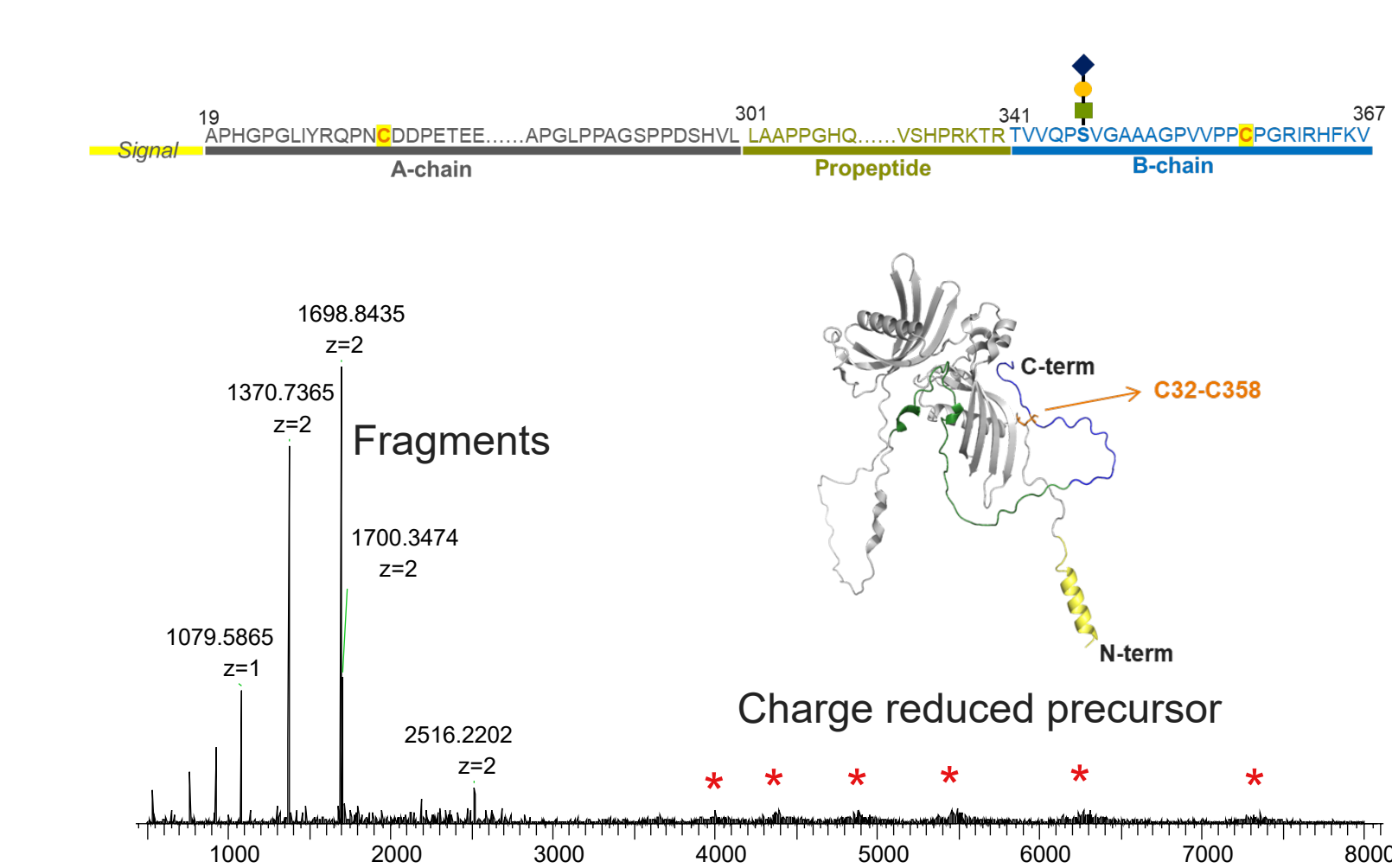
Figure 6. Direct Mass Technology mode Analysis of SARS-CoV-2 Spike Variants



## Native top-down

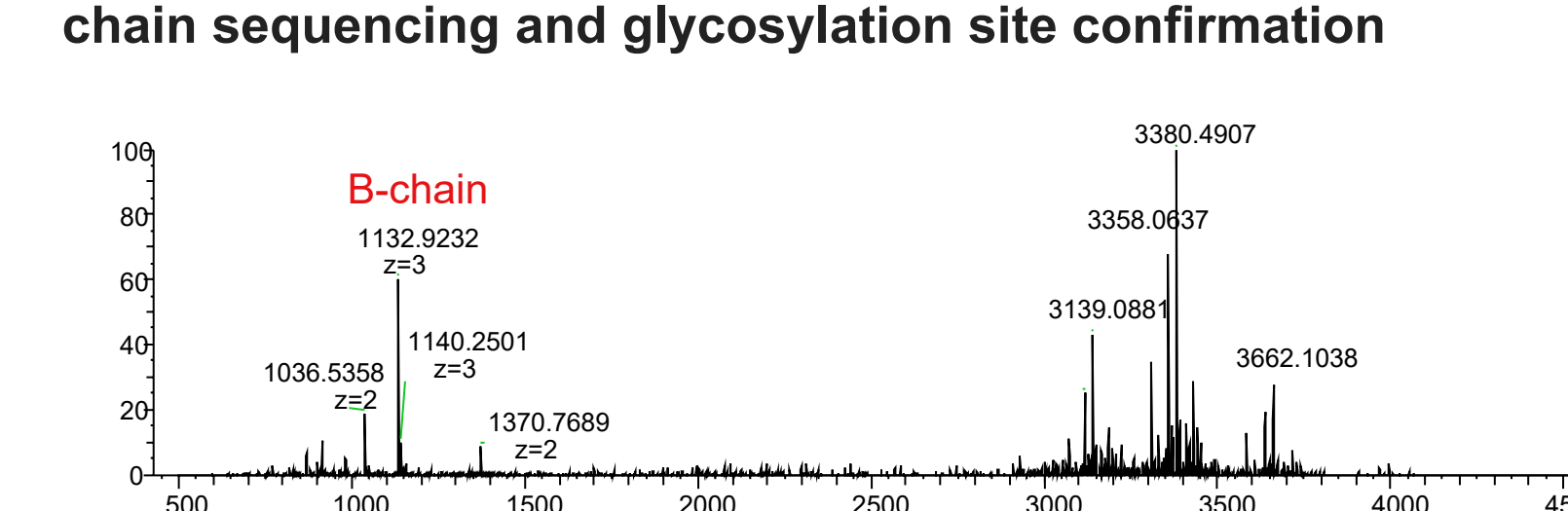
- Native top-down analyses using Hybrid fragmentation confirm sequence, resolve chain connection, and identify glycosylation site

Figure 7. EThcD spectrum and sequence map to resolve chain connection of Fetuin A



Although N terminal of B-chain is connected to the C-terminus of A-chain through a propeptide, EThcD fragments demonstrate the B-chain is disulfide bonded to the N-terminal of A-chain through cys32-cys358

Figure 8. UVPD spectrum and fragments map for Fetuin B-chain sequencing and glycosylation site confirmation



- UVPD top-down analysis could provide 95% sequence and identify O-glycosylated site at Ser346.

## Conclusions

- By combining different MS methods such as I2MS, PTCR and EThcD, we could successfully characterize heavily glycosylated Fetuin.
- Native MS combined with PTCR could resolve proteoforms overlapping in m/z to identify more proteoforms.
- Top-down analyses using hybrid fragmentation confirmed the sequence and PTM site.
- Direct Mass Technology not only isotopically resolved numerous proteoforms of Fetuin but also determine charge states, thus MW of Spike proteins

## References

1. Huguet R et al, Anal.Chem.2019,15732-15739
2. Lin, Yu-Hsien, Vojtech Franc, and Albert JR Heck. Journal of proteome research 17.8 (2018): 2861-2869

## Acknowledgements

1. Christopher Mullen, Mike Senko, Kristina Szreni?c, Daniel Hermanson, Kyle Fort, Florian Grosse-Goosmann from Thermo Fisher Scientific
  2. Ryan Feller, Ken Durbin from Proteinaceous
  3. Prof. Ganesh Anand, Theresa S.C. Buckley, Sean M. Braet from Penn State University
- Jared Kafader, Prof. Neil Kelleher from Northwestern University

## Trademarks/licensing

© 2022 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.

- Direct Mass Technology mode provides the charge states and calculation of the molecular weight profile of each variant
- Multiple forms of HSP70, 90 and cell matrix proteins were detected
- Multiple confirmations of Spike trimers were observed