

ECD and EID fragmentation of peptides and intact proteins using a Quadrupole Time-of-Flight Mass Spectrometer

Valery Voinov^{1,2}, Yury Vasilev^{1,2}, Nathan Lopez^{1,2}, Joseph Beckman¹, Michael Knierman³, Christian Klein⁴, Kenneth Newton⁴, Ruwan Kurulugama⁴, George Stafford⁴, and John Fjeldsted⁴

¹ LPI, Oregon State University, Corvallis, OR ² e-Msion Inc., Corvallis, OR ³ Eli Lilly and Company, Indianapolis, IN ⁴ Agilent Technologies, Santa Clara, CA

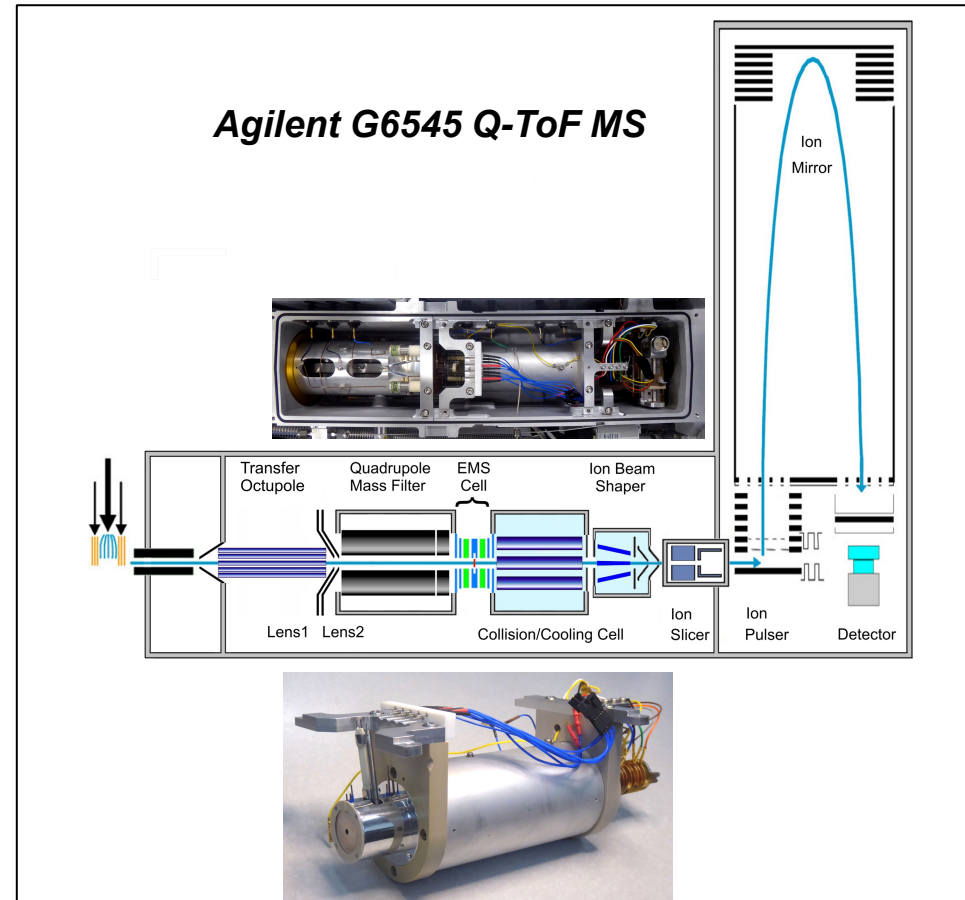


Figure 1. ECD cell and its position in 6545 Q-ToF

- Fort et al. **WOH pm 02:50** "Implementing an Electrostatic ECD Cell on a Q-Exactive Enabling ECD and EChcD Fragmentation" *ASMS 2018 San Diego, CA, June 03-07.*
- Vasil'ev et al. **ThP 800** "Hybrid ECD Methods for Middle-Down And Top-Down Proteomics Implemented in a Benchtop Quadrupole Orbitrap Mass Spectrometer" *ASMS 2018 San Diego, CA, June 03-07.*
- Hill et al. **TOB am 09:30** "On-Line Nanolc-Ion Mobility-Electron Capture Dissociation Tandem MS Analysis of Peptide Mixtures and Glycoprotein Digests on an IMQTOF Mass Spectrometer" *ASMS 2018 San Diego, CA, June 03-07.*

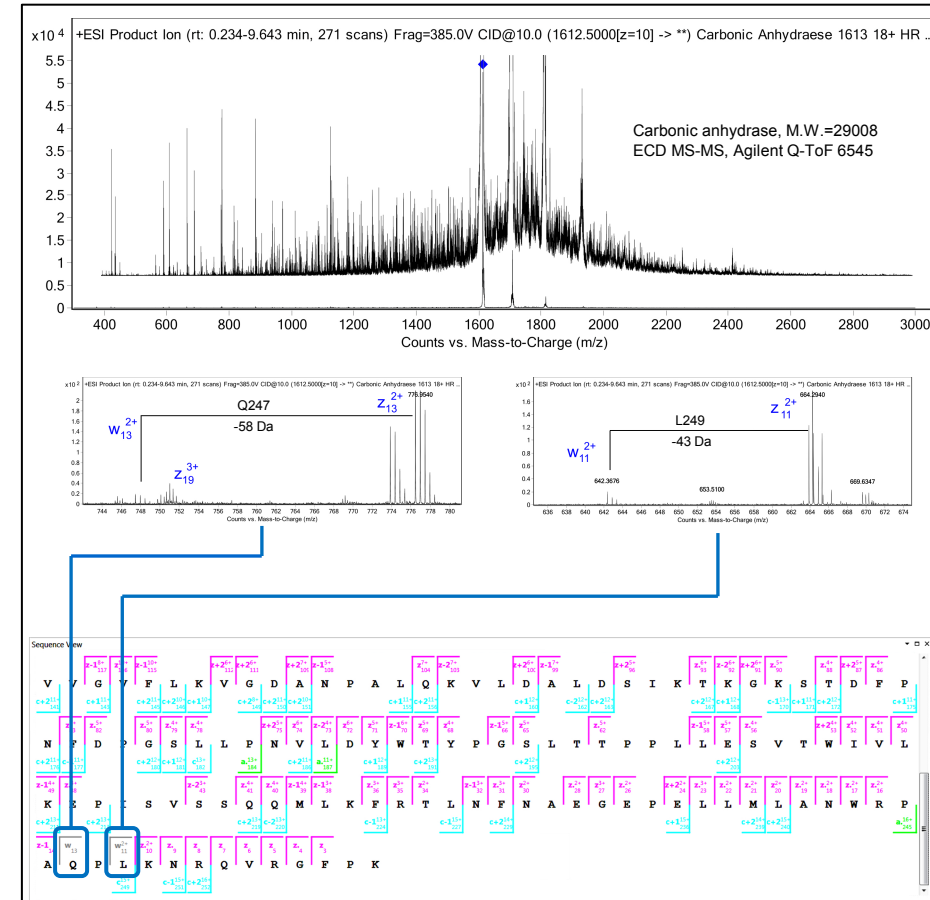


Figure 2. ECD MS-MS spectrum and sequence coverage and fragment matching at 5 ppm mass error for Carbonic Anhydrase 18+ precursor. Insert shows w ions.

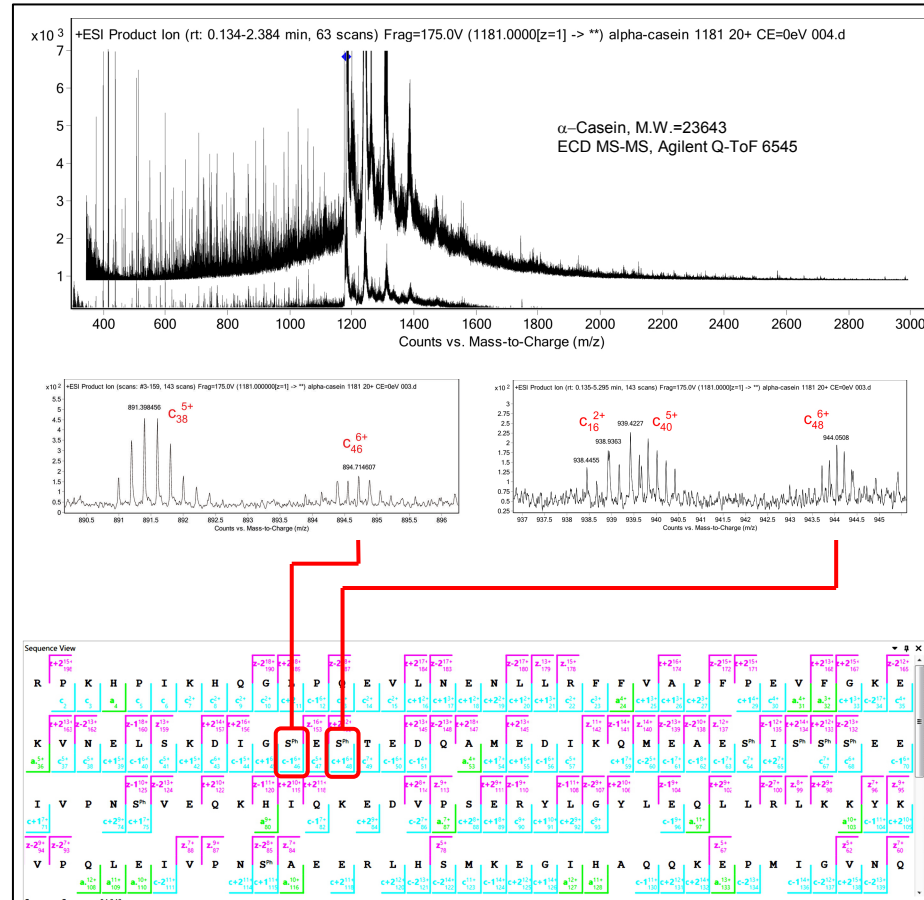


Figure 3. ECD MS-MS spectrum and sequence coverage at 5 ppm mass error for α -casein 20+ precursor. Insert shows c-ions defining phosphorylation sites.

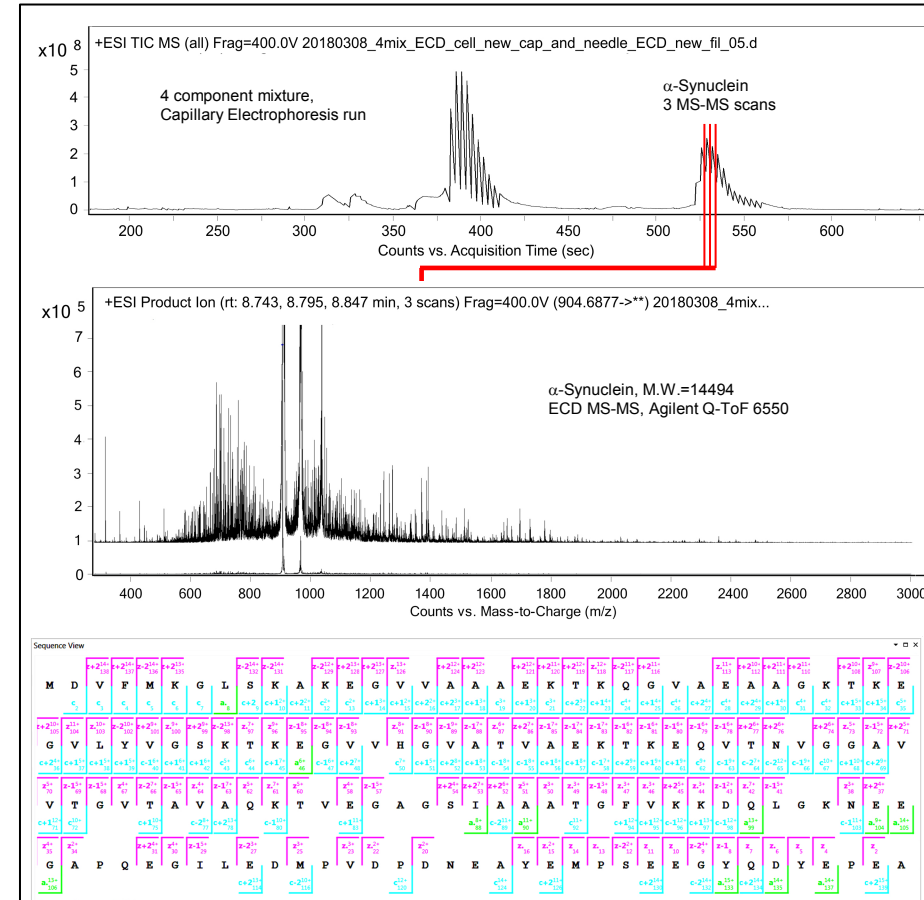


Figure 4. Capillary Electrophoresis chromatogram, ECD MS-MS spectrum and sequence coverage and fragment matching at 10 ppm mass error for α -synuclein.

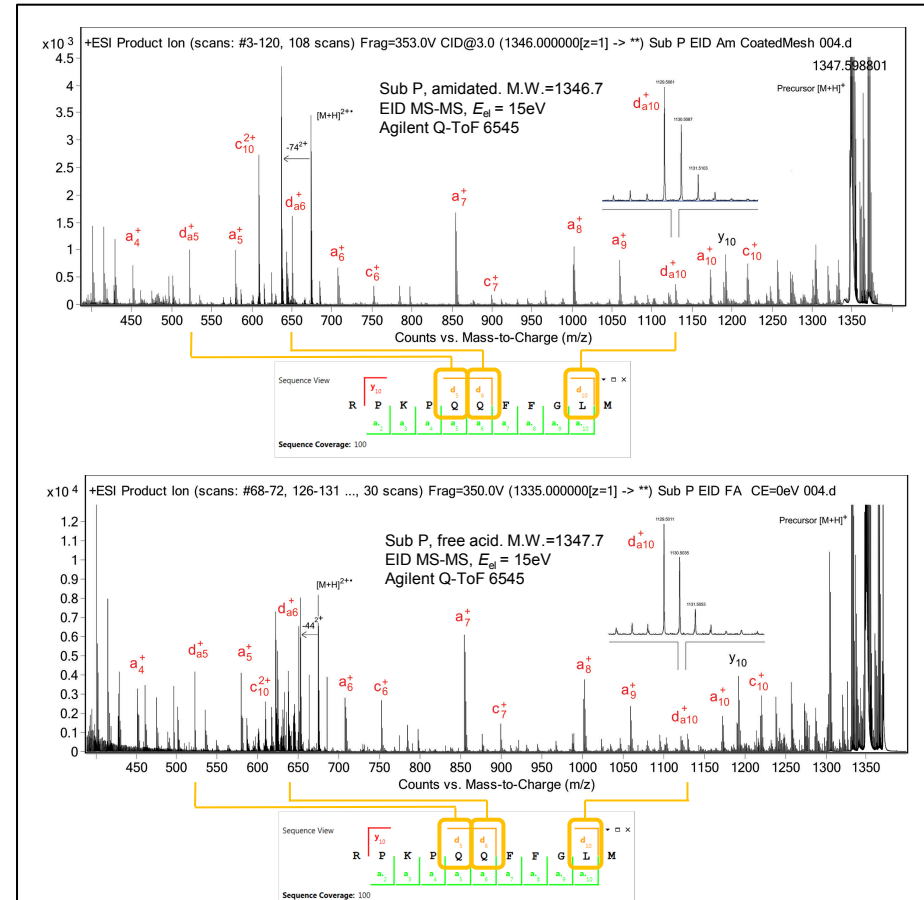


Figure 5. EID (Electron Ionization Dissociation) product ion spectra of Amidated (top) and Free Acid (bottom) forms of substance P. Inserts show d ions.

e-Msion ECD technology:

- Compatible with virtually any mass spectrometry platform (Fig.1);
- Original functionalities of the retrofitted instruments are retained;
- Allows Leu & Ile to be distinguished (Fig.2, 5);
- Preserves Post Translational Modifications (Fig.3);
- Fast, compatible with HPLC, CE and Ion Mobility (Fig.4);
- Higher electron energy modes available (Fig.5);
- Provides high sequence coverage;
- Provides low false discovery rate.

Acknowledgement

V.G.V., Y.V.V., N.I.L. and J.S.B. are cofounders of e-Msion, Inc., which is based on technology developed and licensed from Oregon State University. As employees of Oregon State University, a potential conflict of interest exists. This work was supported by the NIH (SBIR 123855). The authors thank Larry M. Nelson (OSU, Corvallis, OR) for fabricating components of the EMS ECD cell. LCMS Spectator (Park, J. et al. *Nat. Methods* 2017, 14, 909–914) was used for data analysis.

