

# A Fast and Efficient Arginine-Specific Protease for Proteomic Applications

Chris Hosfield<sup>1</sup>, Jessica Wohlfahrt<sup>2</sup>, Jennifer Guergues<sup>2</sup>, Alba Katiria Gonzalez Rivera<sup>1</sup>, Ethan Strauss<sup>1</sup>, Michael Rosenblatt<sup>1</sup>, Marjeta Urh<sup>1</sup> & Stanley Stevens Jr.<sup>2</sup>

<sup>1</sup>Promega Corporation, 2800 Woods Hollow Rd, Madison, WI 53711; <sup>2</sup>University of South Florida, 4202 E Fowler Ave, Tampa, FL, 33620

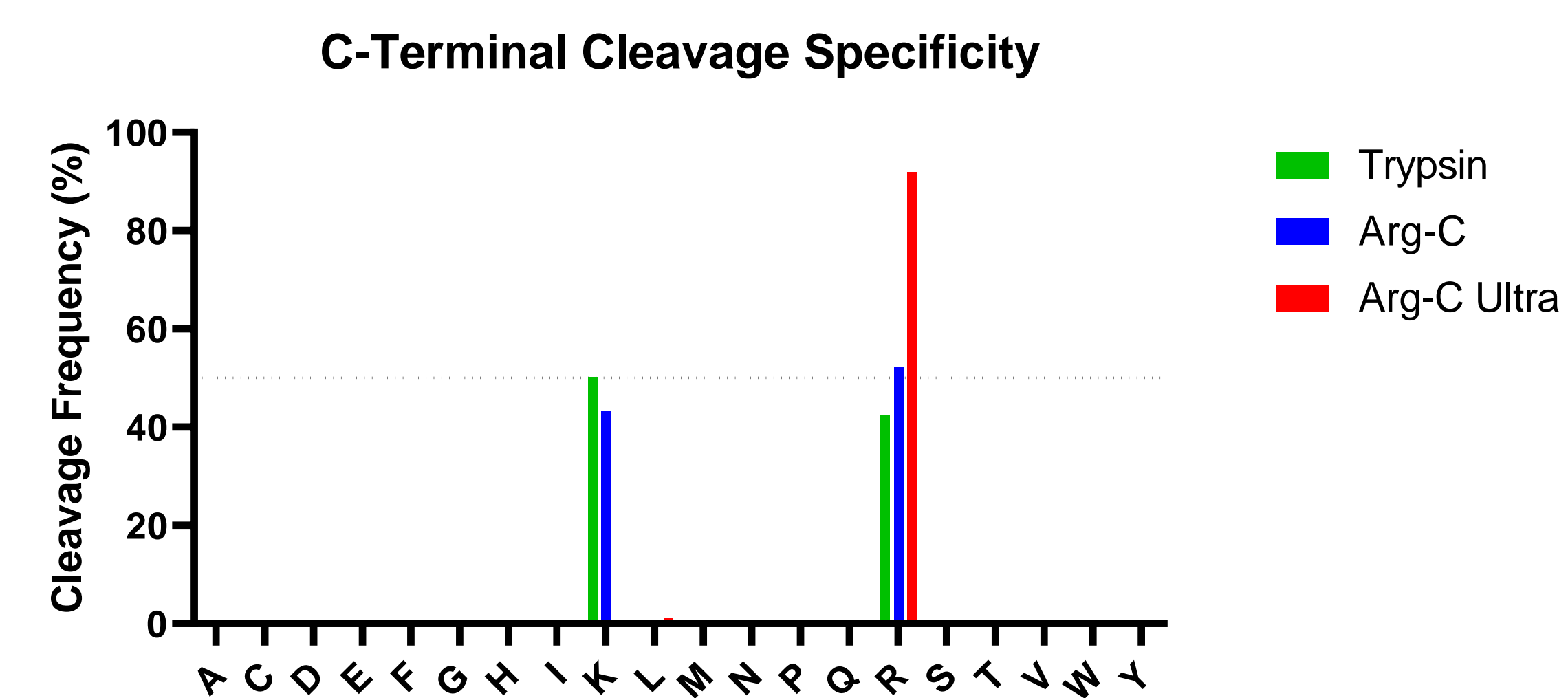


## 1. Introduction

In bottom-up proteomic studies, poor digestion due to low specificity and efficiency complicates LC-MS/MS analysis by increasing sample complexity and reducing sensitivity. Aside from Trypsin and Lys-C, most commonly used proteases suffer from either poor efficiency or specificity.

Here we utilize an arginine-specific protease which is both highly efficient and specific, unlike the commonly used Arg-C protease which cleaves extensively at lysine residues. The improved protease, called Arg-C Ultra, was successfully used in a variety of applications including characterization of human cell extracts, antibodies and post-translational modifications of histones.

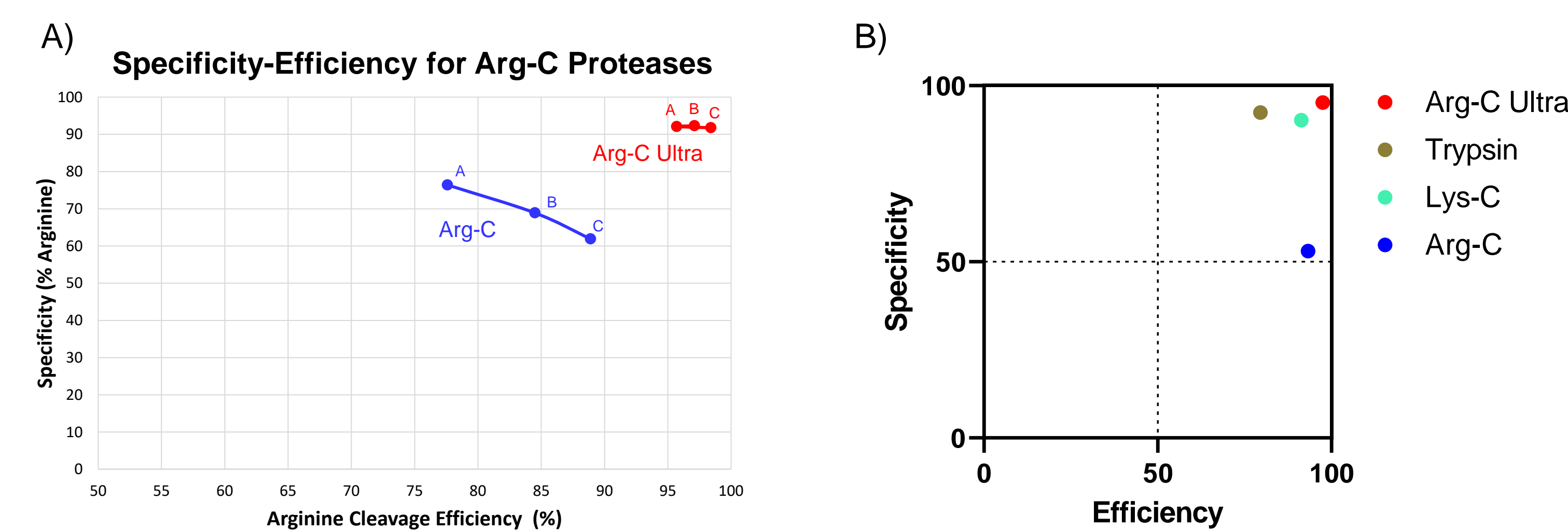
## 2. Arg-C Ultra has Significantly Improved Specificity



• Human K562 cell extract was digested overnight at 1:50 enzyme:substrate ratio.  
• Data were analyzed on an Orbitrap Exploris 240 and searched with Byonic (no enzyme specified)

➤ The digestion specificity of Arg-C Ultra is far superior to standard Arg-C.

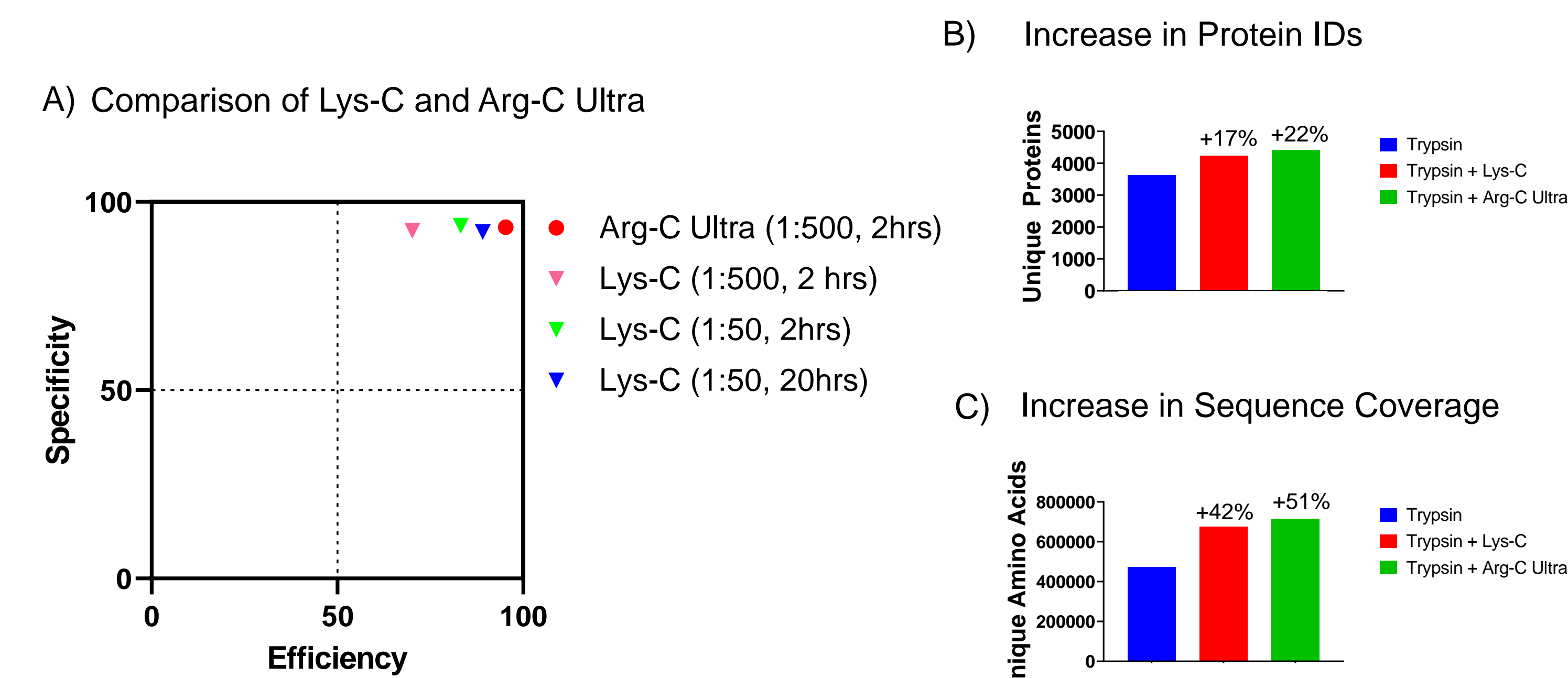
## 3. Arg-C Ultra is a Fast and Efficient Protease



A) Human K562 extract was digested with Arg-C or Arg-C Ultra for 2 hours at 1:500 (A), 1:150 (B) and 1:50 (C). B) K562 extract was digested 1:50 overnight with various proteases.

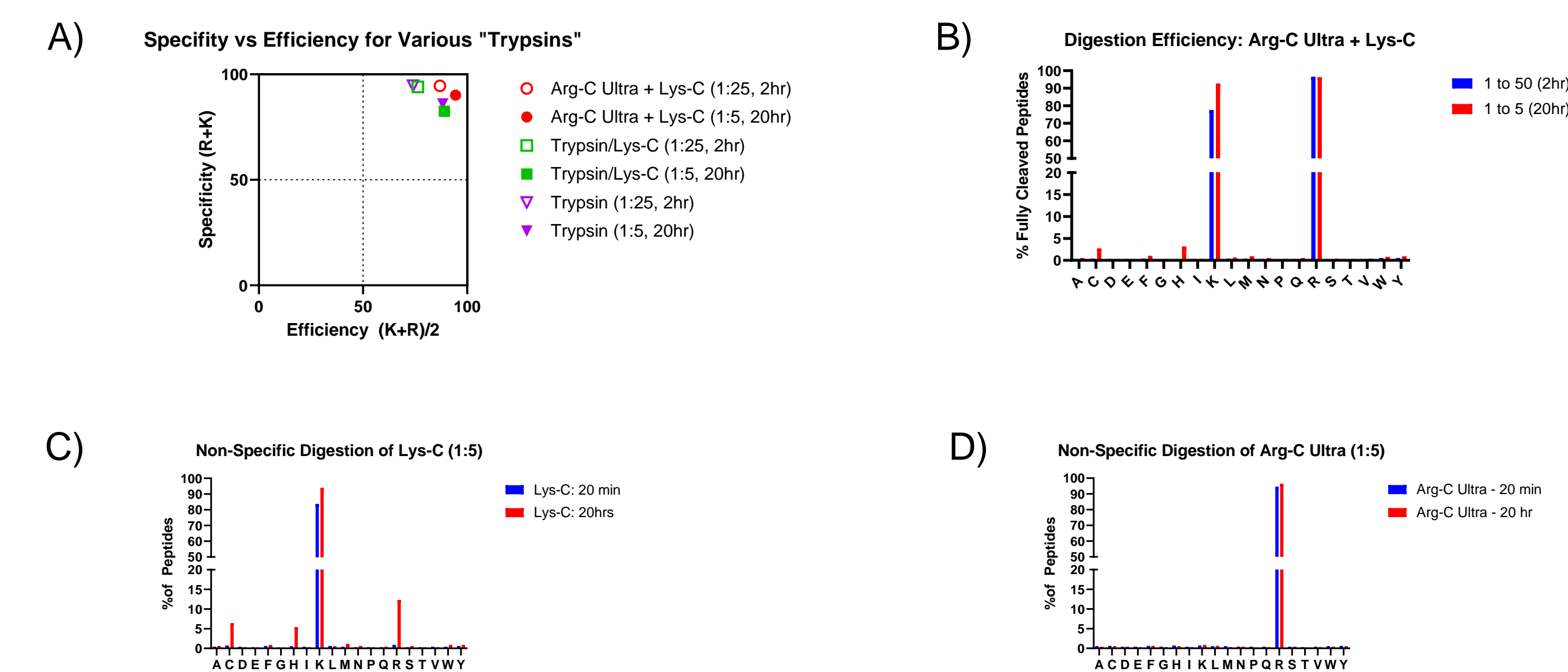
➤ Arg-C Ultra has exceptional digestion efficiency even with short digests using little enzyme.  
➤ Arg-C Ultra has the best combination of efficiency and specificity of MS-grade proteases.

## 4. Arg-C Ultra Outperforms Lys-C



➤ Arg-C Ultra achieves higher digestion efficiency than Lys-C, even with 10X less enzyme and digestion time.  
➤ Arg-C Ultra improves protein IDs and sequence coverage relative to Trypsin to a greater extent than Lys-C.

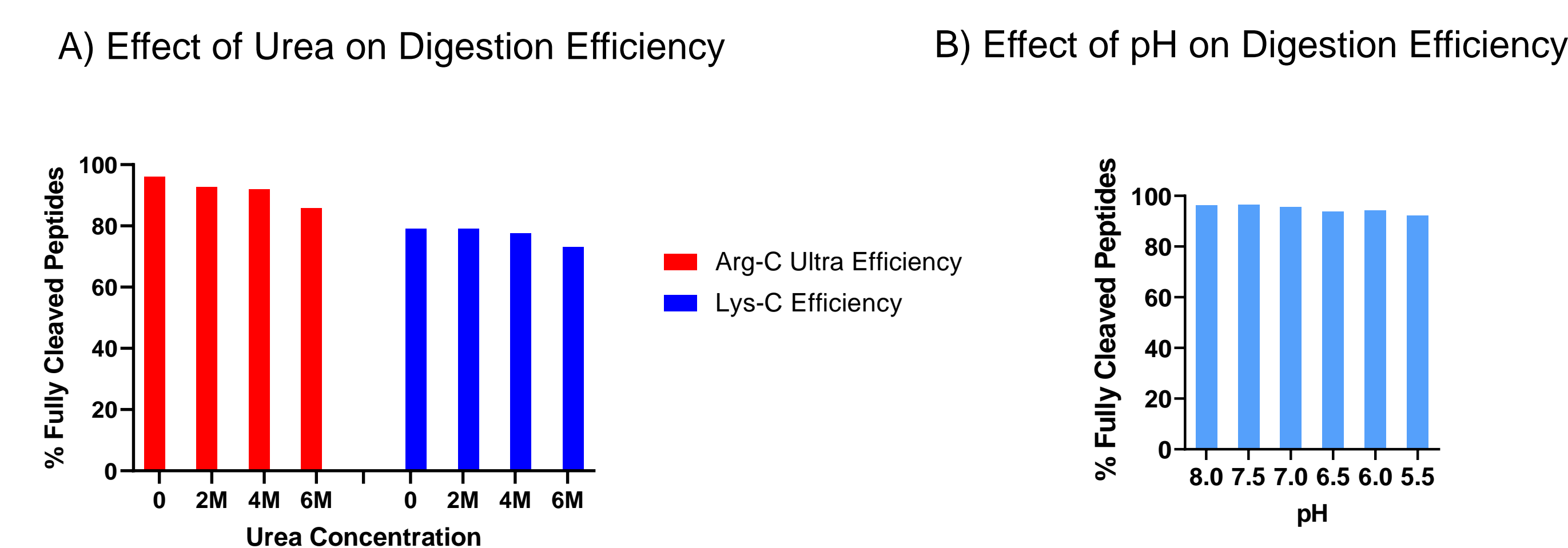
## 5. Arg-C Ultra + Lys-C = "Trypsin Ultra"



Human K562 extract was digested with either trypsin, Trypsin/Lys-C or a mixture of Lys-C and Arg-C Ultra. Data were searched with Byonic (No enzyme specified)

➤ Combining Arg-C Ultra with Lys-C results in better performance than with Trypsin or Trypsin/Lys-C.  
➤ Arg-C has higher digestion specificity than Lys-C.

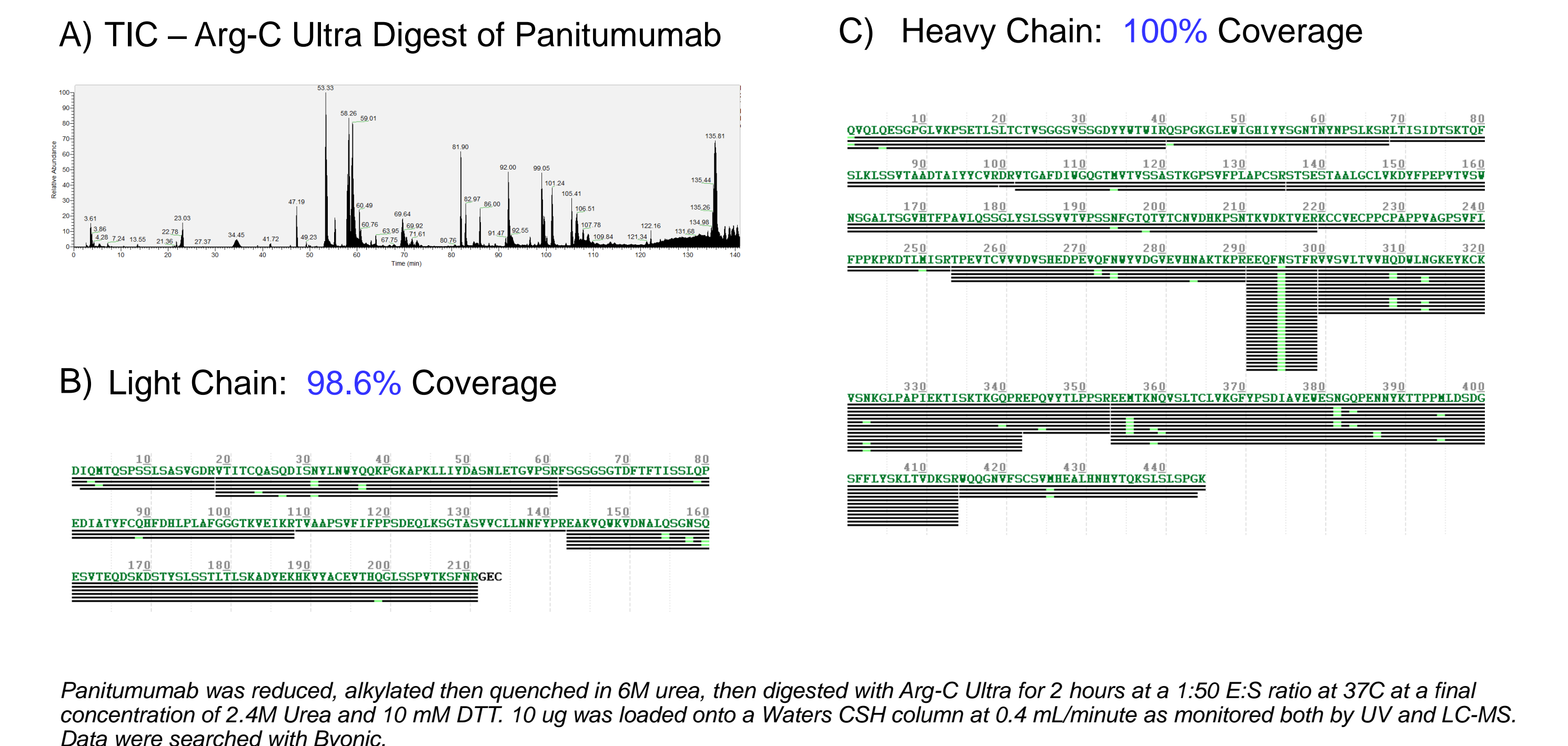
## 6. High Digestion Efficiency both at Low pH or High Urea



Human K562 extract was digested with Arg-C or Lys-C at 1:100 for 2 hours at 37C at various urea concentrations or pH values. Data were searched with Byonic (No enzyme specified)

➤ Arg-C Ultra is more efficient in 6M urea than Lys-C.  
➤ Arg-C Ultra is highly efficient over a wide pH range (~5-9).

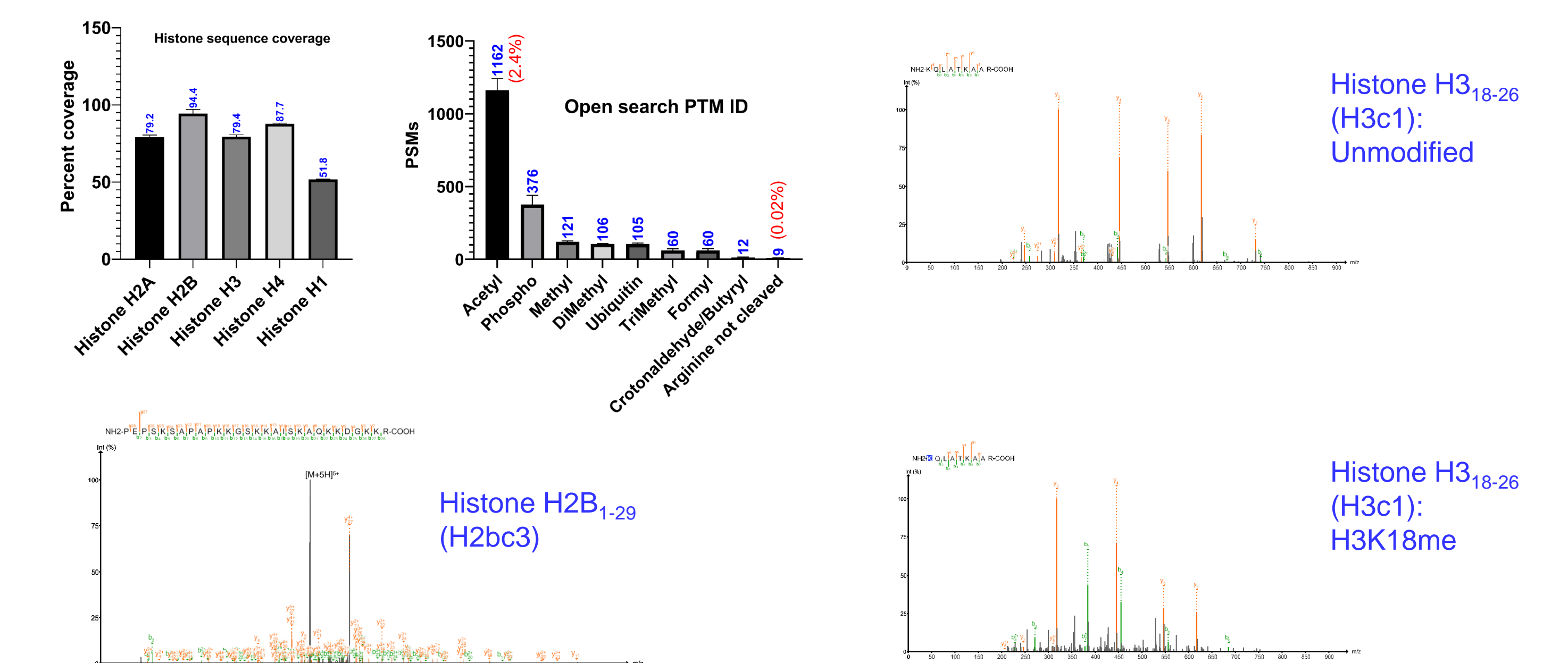
## 7. Highly Specific Peptide Mapping of Therapeutic IgGs



Panitumumab was reduced, alkylated then quenched in 6M urea, then digested with Arg-C Ultra for 2 hours at a 1:50 E:S ratio at 37C at a final concentration of 2.4M Urea and 10 mM DTT. 10 ug was loaded onto a Waters CSH column at 0.4 mL/minute as monitored both by UV and LC-MS. Data were searched with Byonic.

➤ High sequence coverage of Panitumumab obtained after digestion with Arg-C Ultra for 2 hours.

## 8. Histone Analysis: High Sequence Coverage and PTM ID



Histones were extracted from mouse microglia with the EpiQuik Total Histone Extraction Kit, digested with ArgC-Ultra, and analyzed via DDA-PASEF on a timsTOF Pro. MS Fragger search results show high sequence coverage (spanning the N-terminal tail) and PTM ID.

➤ ArgC-Ultra digestion allows detection of PTM-containing and unmodified cognate histone peptides for accurate histone PTM quantitation.

## 9. Advantages of Arg-C Ultra

- Highest performing MS-grade protease
  - Specific for arginine.
  - Efficient: < 5% missed cleavages even with 1:500, 2hr digest.
- When combined with Lys-C, Arg-C Ultra + Lys-C is superior to trypsin.
- Superior to Lys-C:
  - Higher digestion efficiency despite requiring less enzyme.
  - Higher specificity.
  - More orthogonal to Trypsin for protein coverage in K562 digest.
  - More efficient even in 6M urea.
- Wide pH range, potentially useful for peptide mapping at low pH.
- Useful for producing high coverage maps of therapeutic proteins.
- Useful for comprehensive PTM mapping as evidenced by Histone analysis.
  - Particularly helpful for monitoring PTMs on or around Lysine residues.