

Enhancing trypsin digestion with Lys-C and Arg-C proteases

Sergei V. Saveliev¹, Laurie Engel¹, Mike Rosenblatt¹, Richard Jones², Michael Ford², Ravi Amunugama², Dave Allen² and Marjeta Urh¹

¹Promega Corporation, Madison, WI; ²MS BioWorks, LLC, Ann Arbor, MN

email: sergei.saveliev@promega.com

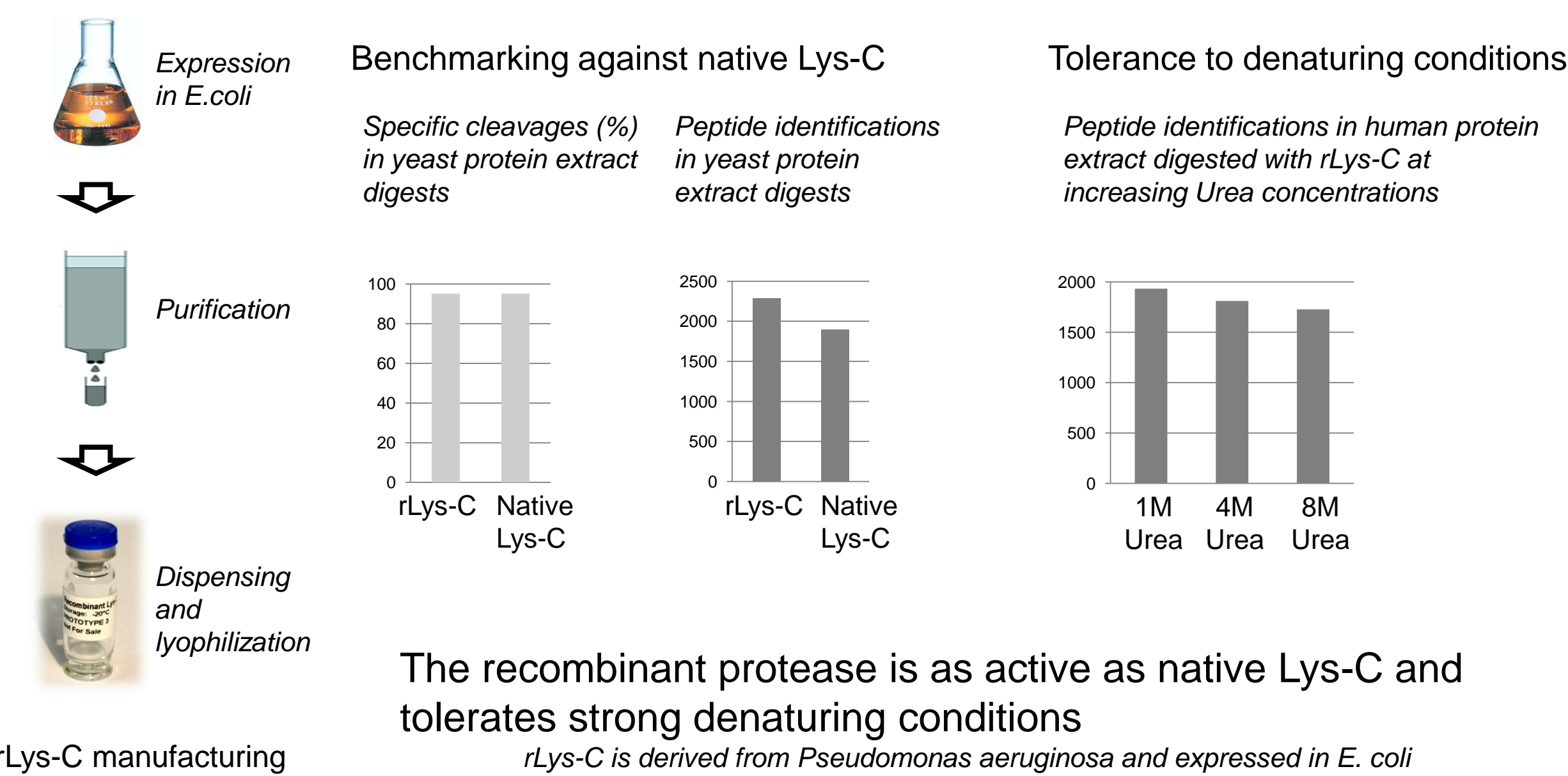


1. Introduction

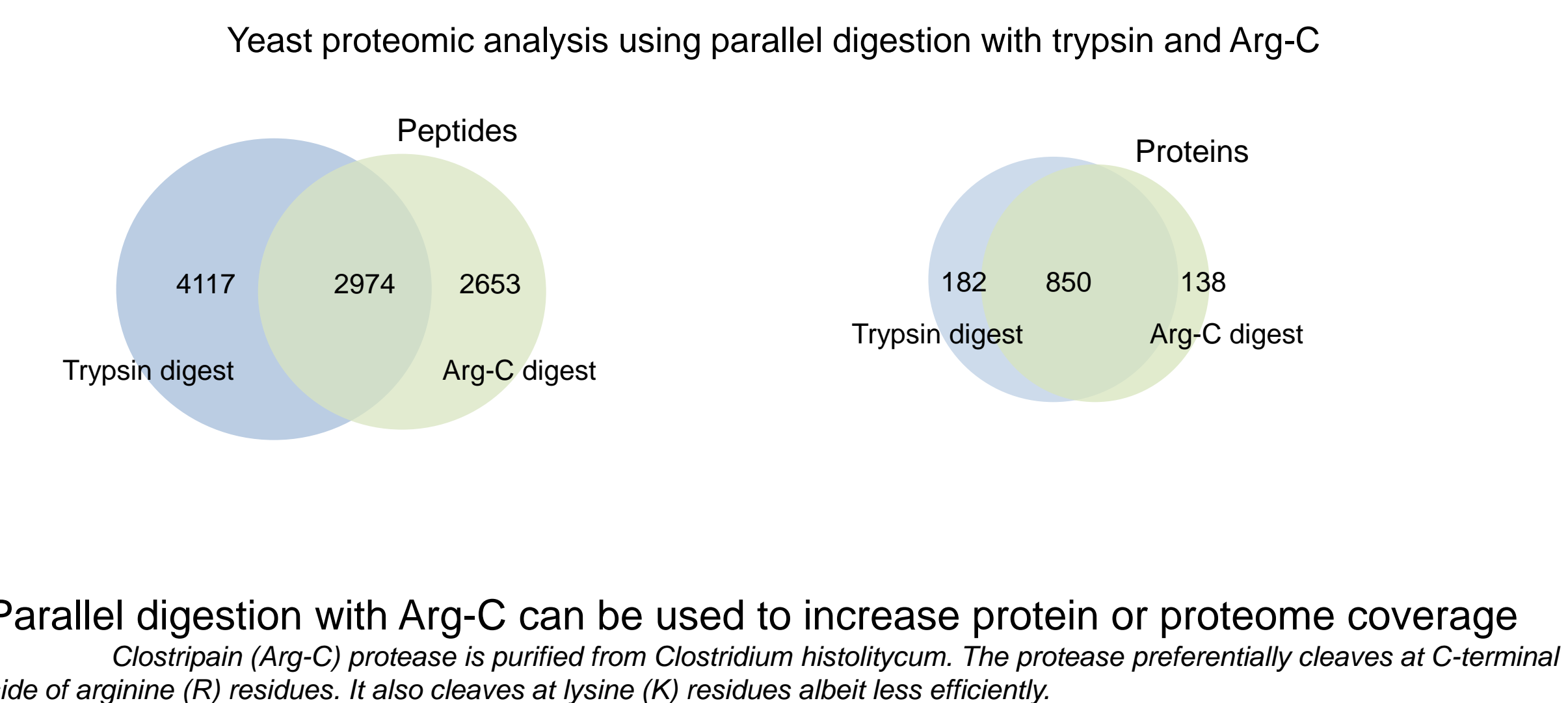
Owing to efficient proteolysis and particular advantages of trypsin-generated peptides, trypsin is the most widely used protease in mass spectrometry. However, it is often overlooked that trypsin has shortcomings such as incomplete proteolysis, poor digestion of tightly folded proteins or an inability to cleave arginine and lysine residues adjacent to proline. These shortcomings negatively affect mass spec analysis, particularly advanced applications such as protein quantitation.

We show here that Lys-C and Arg-C proteases address trypsin shortcomings. By supplementing trypsin with Lys-C we have achieved $\geq 95\%$ digestion efficiency as judged by the level of completely digested peptides. In addition, by utilizing the remarkable resistance of Lys-C to denaturing conditions we have adapted this trypsin/Lys-C mix to digest a proteolytically resistant protein. Finally, we demonstrate the unique utility of Arg-C protease based on its ability to cleave the residues adjacent to proline.

2. Recombinant Lys-C protease

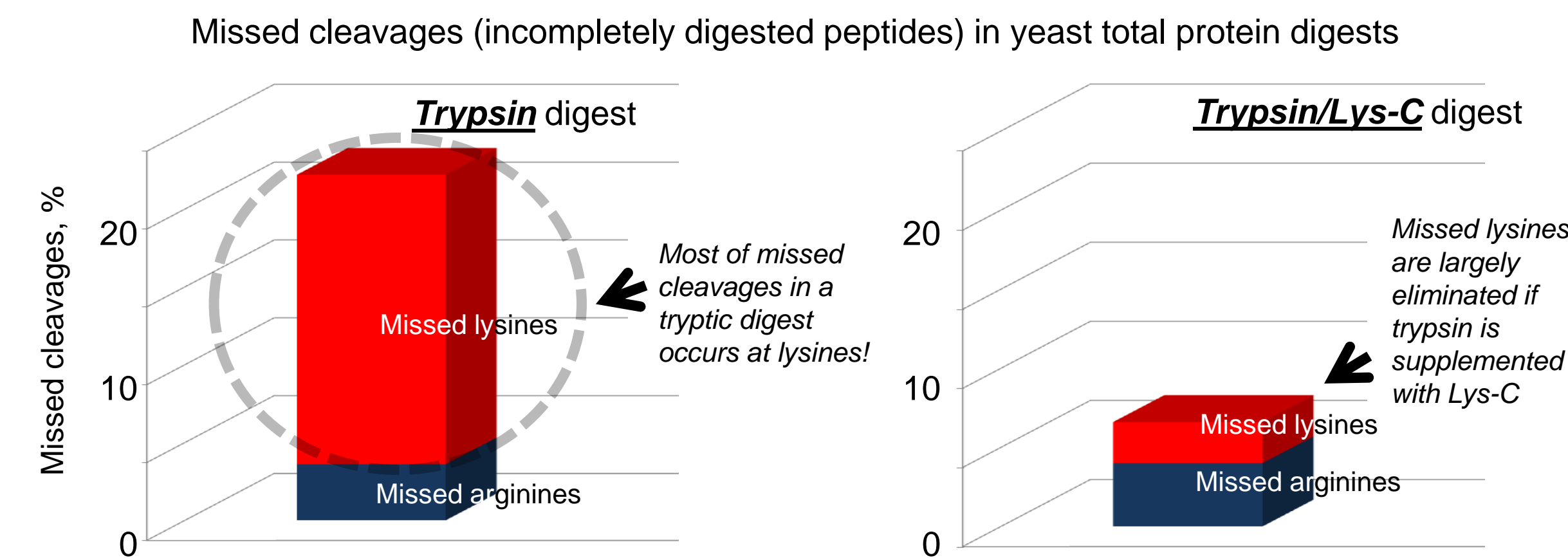


3. Clostripain (Arg-C) protease



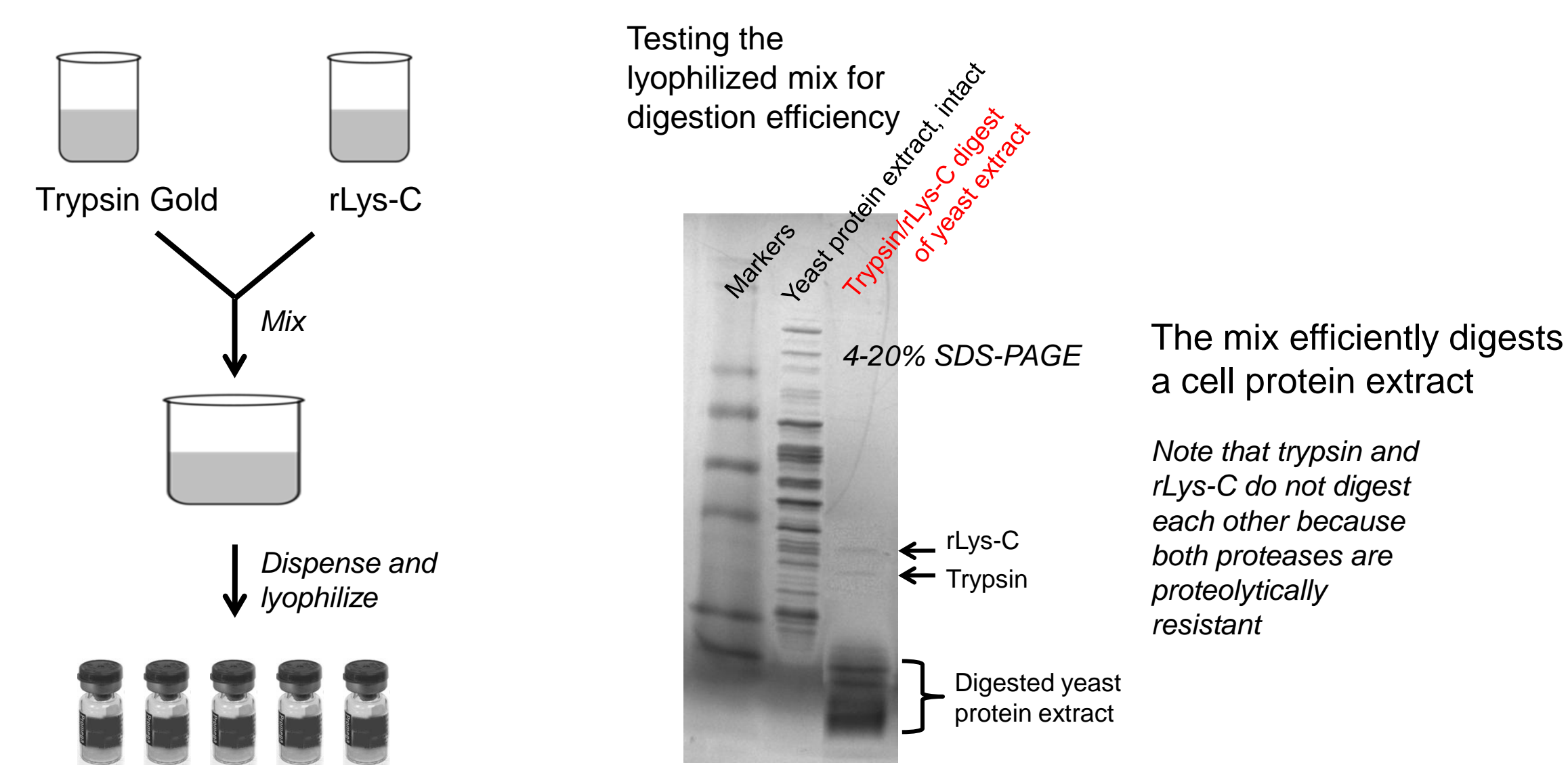
4. Advantage of Lys-C-supplemented trypsin

Typically, ~20-30% peptides in a tryptic digest are incompletely digested. Incomplete digestion is a major source of irreproducibility of mass spec analysis or inaccurate protein quantitation. In contrast to a logical assumption, missed R and K in a tryptic digest are not represented in equal amounts. Most of missed cleavages occur at lysines.



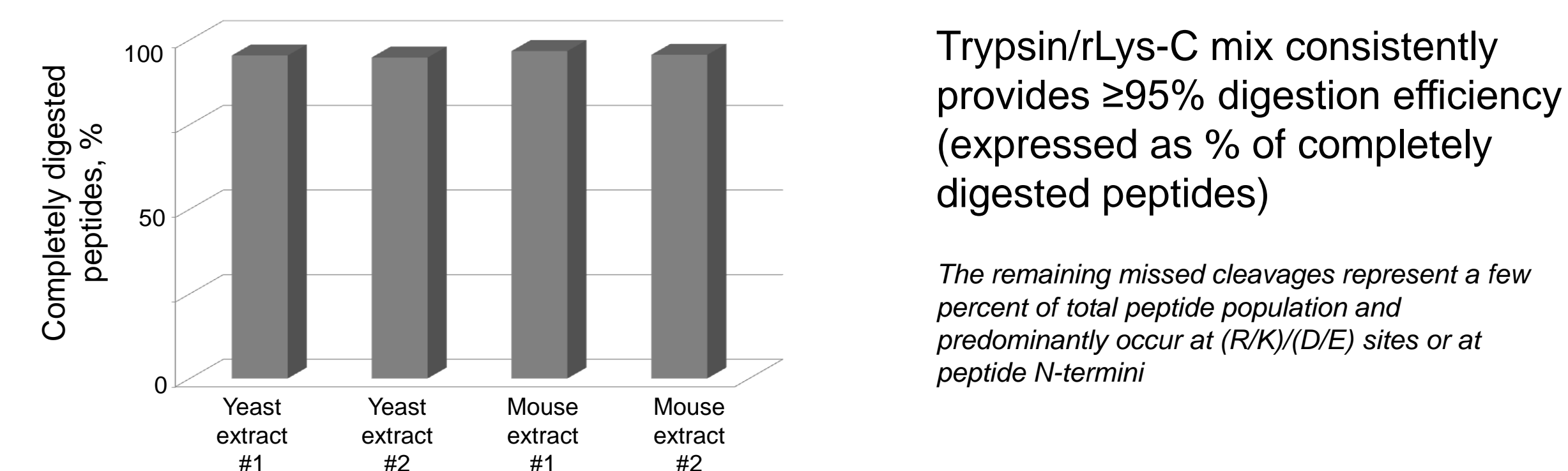
Supplementing trypsin with Lys-C dramatically improves digestion efficiency

5. Preparation of trypsin/rLys-C mix



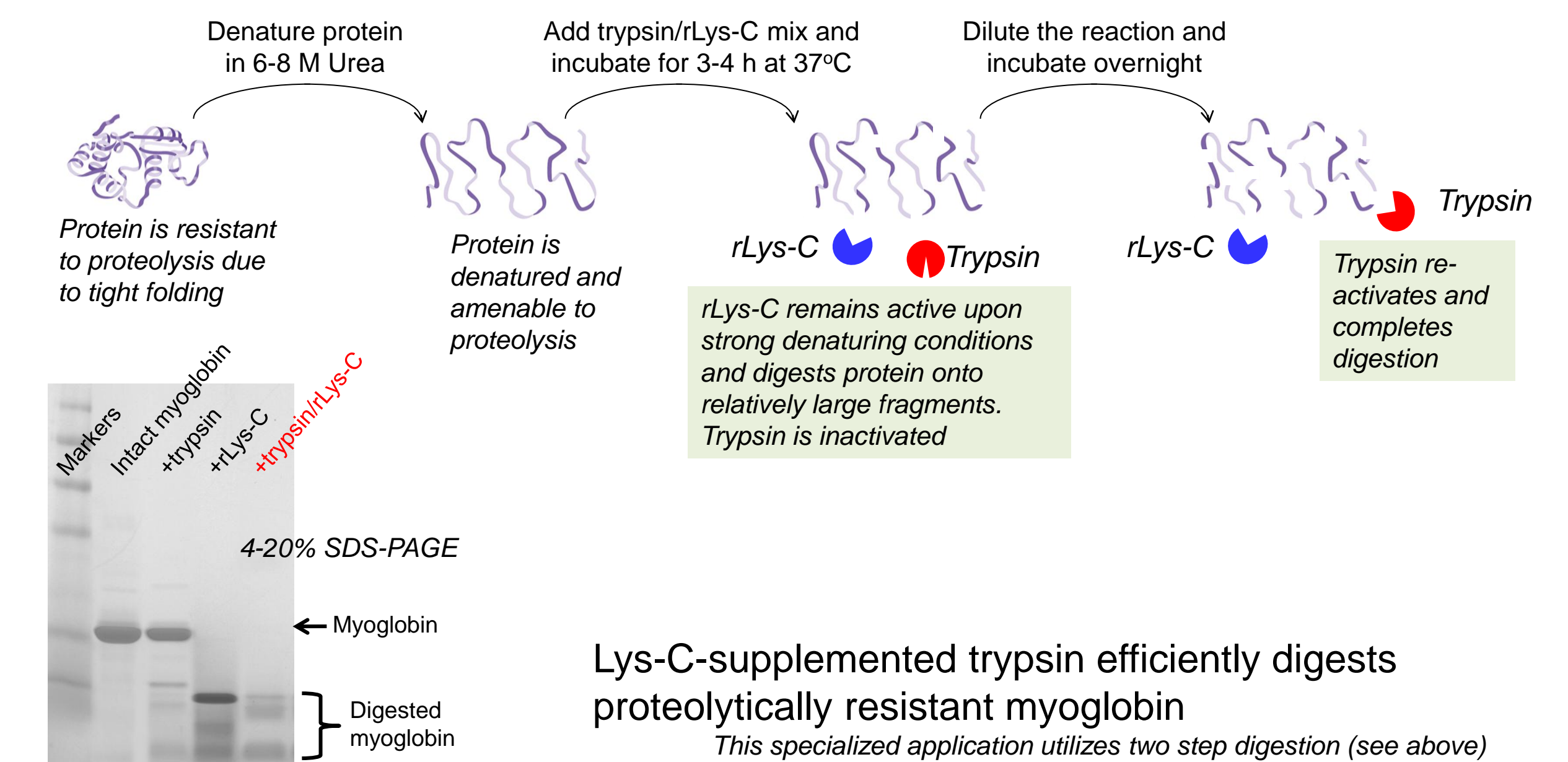
6. Validation of Trypsin/rLys-C mix

Digestion of yeast and mouse protein extracts with Trypsin/rLys-C mix



Experimental conditions: total yeast and mouse protein extracts were digested with trypsin/rLys-C mix at 37°C overnight. The digestion was performed under conventional trypsin-amenable (non- or mildly denaturing) conditions. The digests were analyzed with LTQ Orbitrap Velos

7. Additional, specialized application of Trypsin/rLys-C mix: digestion of tightly folded, proteolytically resistant proteins



8. Arg-C advantage

Cleavage of RP and KP sites with Arg-C in yeast total protein extract

Protease	Unique peptides	Cleavages at arginines	Cleavages at lysines	Cleaved RP sites	Cleaved KP sites
Trypsin	4481	37%*	61%*	0%	0%
Arg-C	3374	56%	41%	97%	44%

*Difference in the share of arginine and lysine cleavages in trypsin digest reflects natural unequal distribution of these residues in proteins rather than trypsin cleavage preference toward lysines

Unlike trypsin, Arg-C efficiently cleaves at arginine and, to a lesser extent, lysine residues adjacent to proline

9. Conclusion

- Supplementing trypsin with rLys-C improves digestion efficiency $\geq 95\%$ peptides are completely digested
- Trypsin/Lys-C mix can also be used to digest tightly folded, proteolytically resistant proteins. This specialized application utilizes Lys-C tolerance to protein denaturing conditions
- Arg-C protease complements trypsin by cleaving arginine and lysine sites before proline. Arg-C can also be used to improve proteome coverage