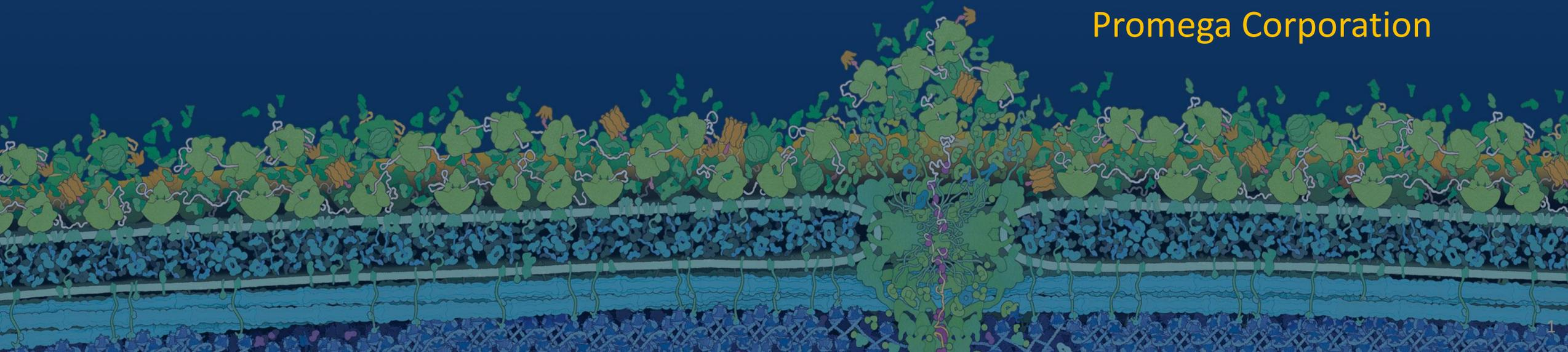
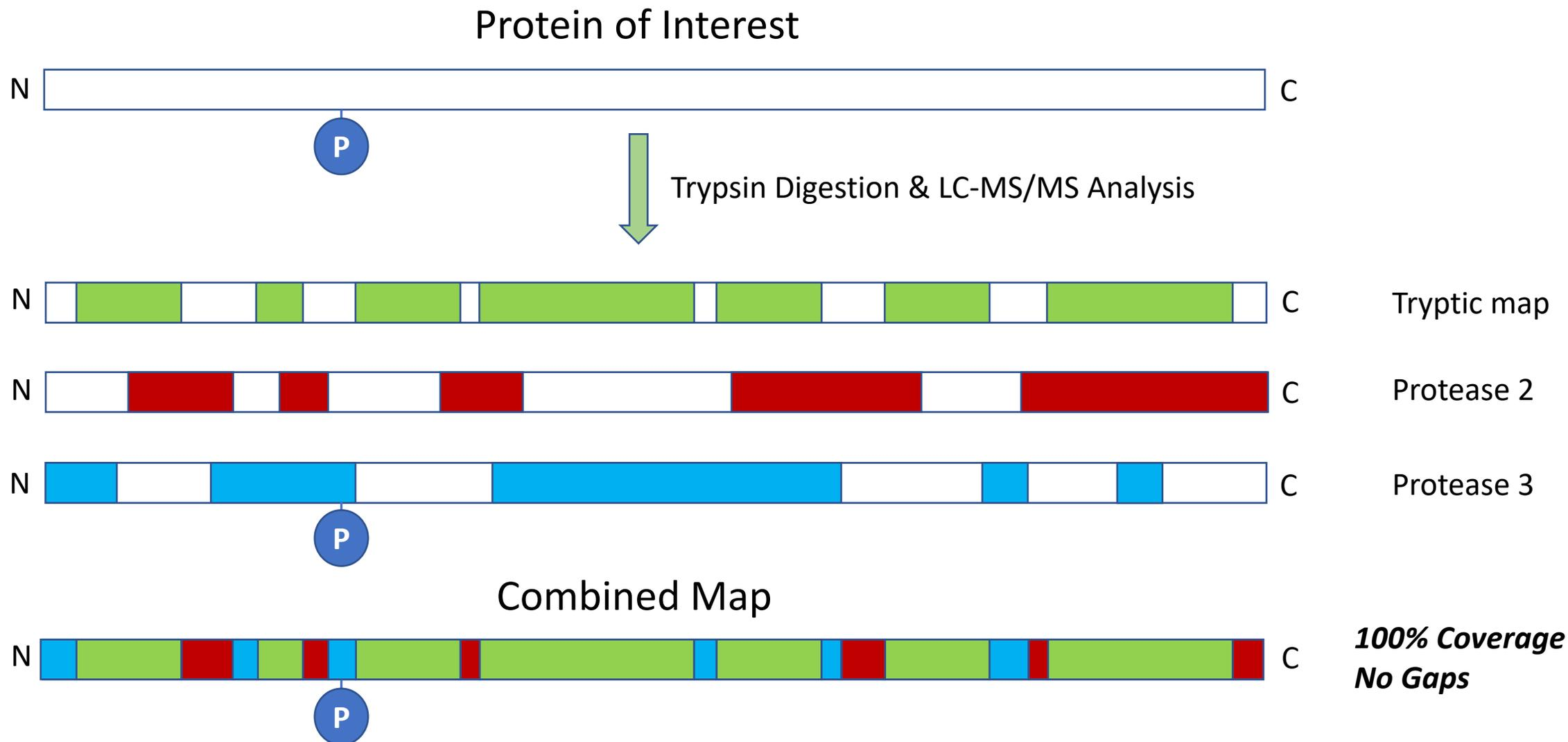


# ProAlanase: A Proline- and Alanine-Specific Protease for Proteomics

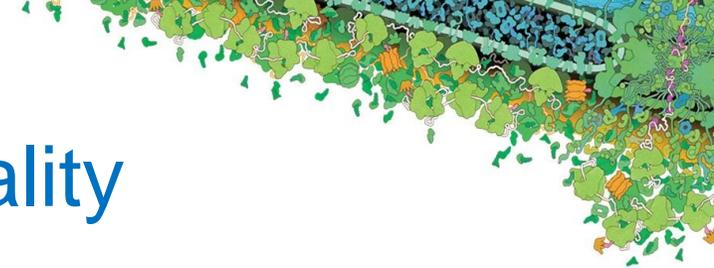
Chris Hosfield, PhD  
Senior Research Scientist  
Promega Corporation



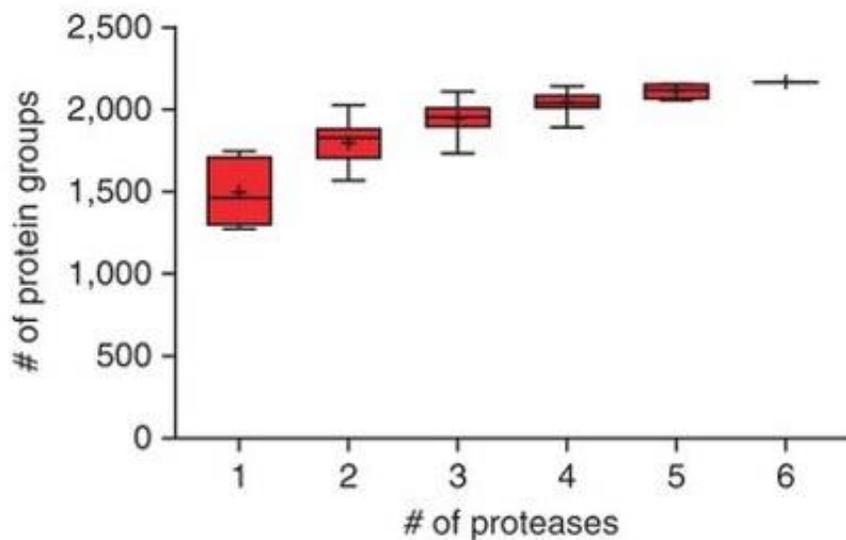
# Trypsin is Great: But Rarely Gives the Full Picture



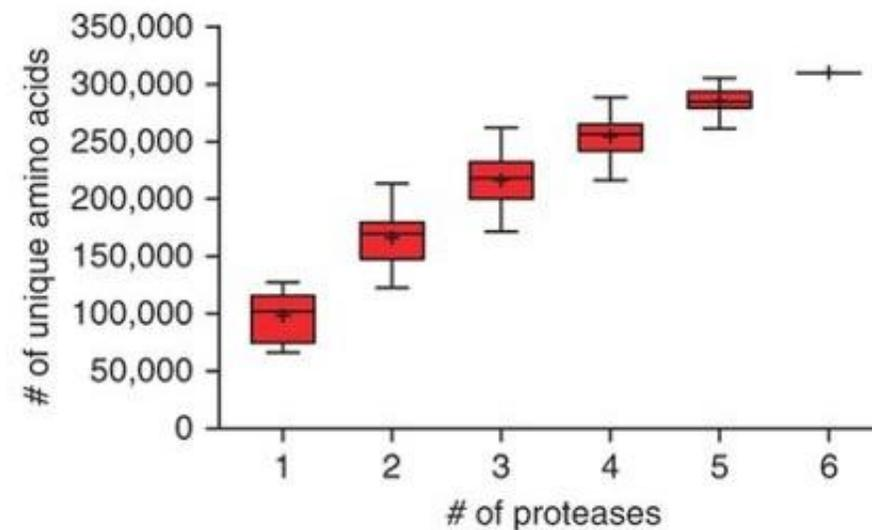
# Using Multiple Proteases Improves Data Quality



Protein Identifications



Amino Acid Coverage



*E. Coli* lysate

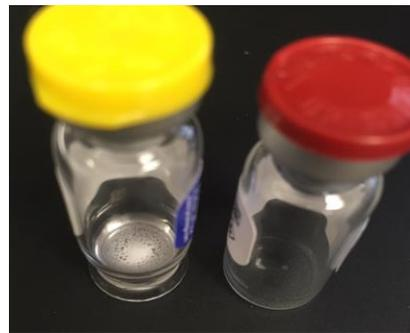
*Giansanti et al. (2016)*  
*Nature Protocols*

# Recent Addition: Recombinant Asp-N (rAsp-N)

- High performance – efficient digestion in one hour
- His-tag for easy removal
- Improved format
  - New formulation enhances product stability and consistency of use
  - V-shaped vial and larger size (10 µg)



V-shape Flat-Bottom

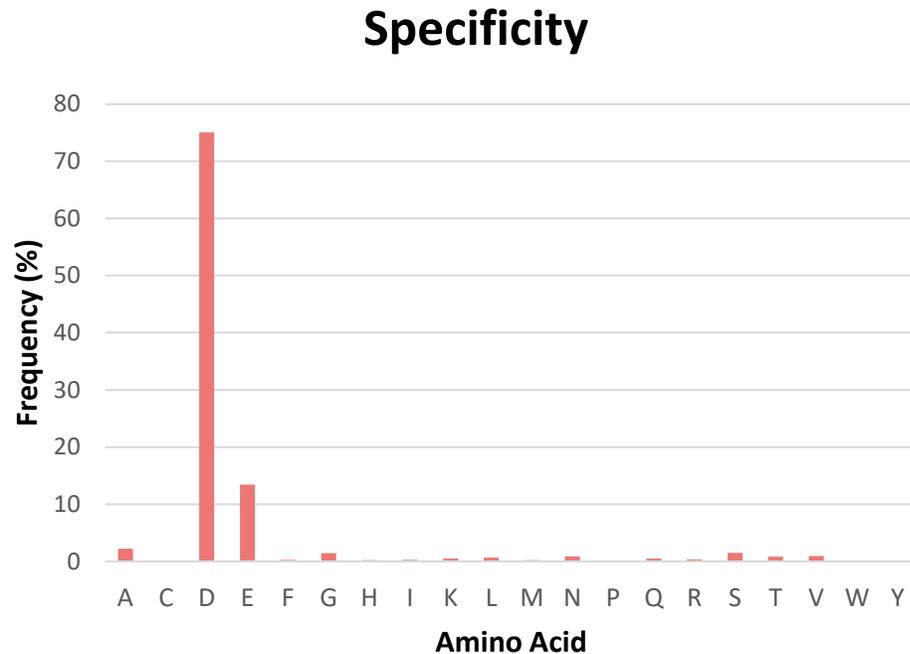


Enhanced formulation

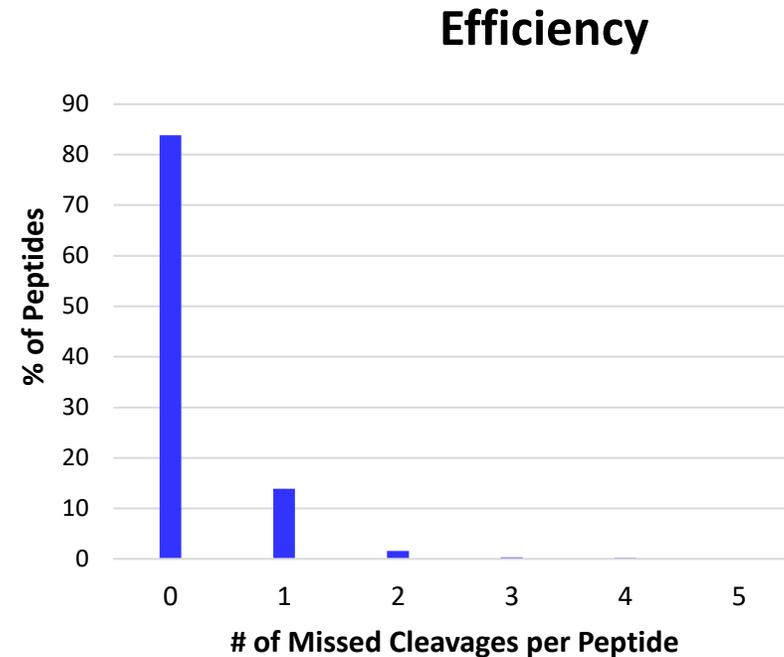
**Better price**  
~1/2 the cost of native per µg

# rAsp-N is a High-Performance Protease

- Yeast extract was digested for **1 hour** at 37°C in ~1.5M urea at a 1:50 ratio

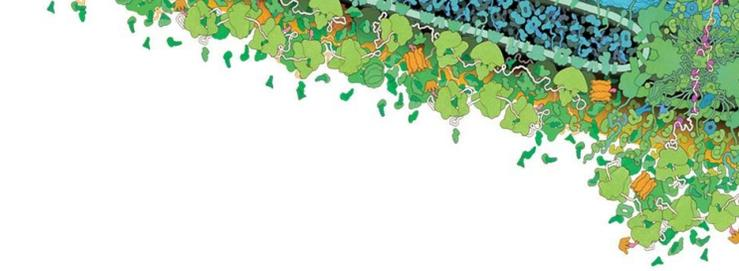


- Primarily cleavage at Asp
- Secondary cleavage at Glu

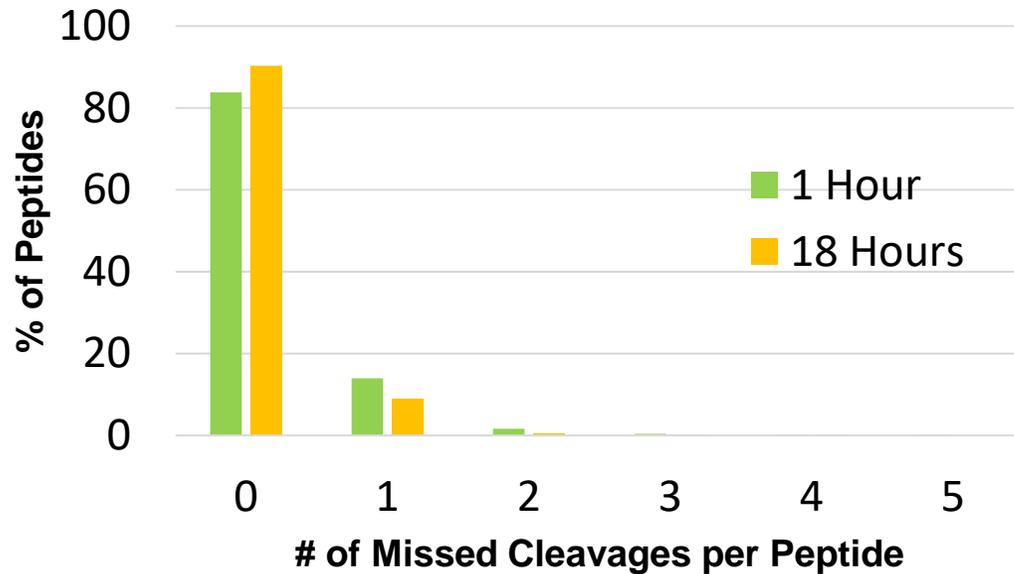


- Cleavage at Asp is ~85% complete

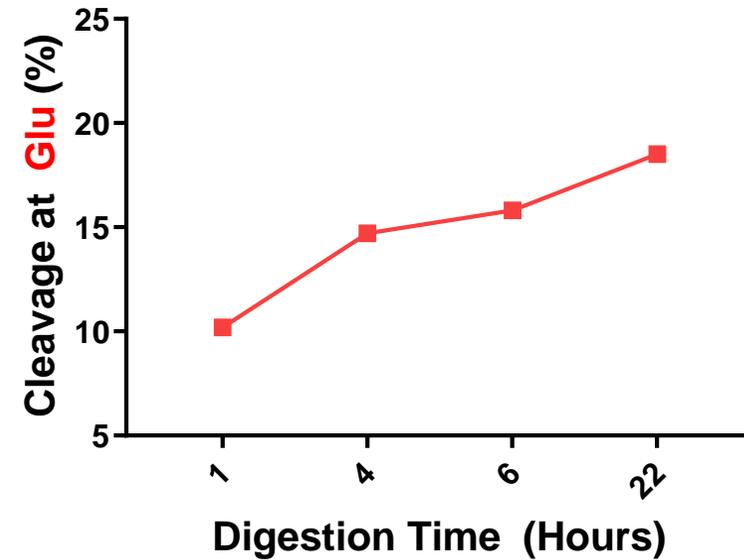
# What Happens in Longer Digests?



### Cleavage Efficiency



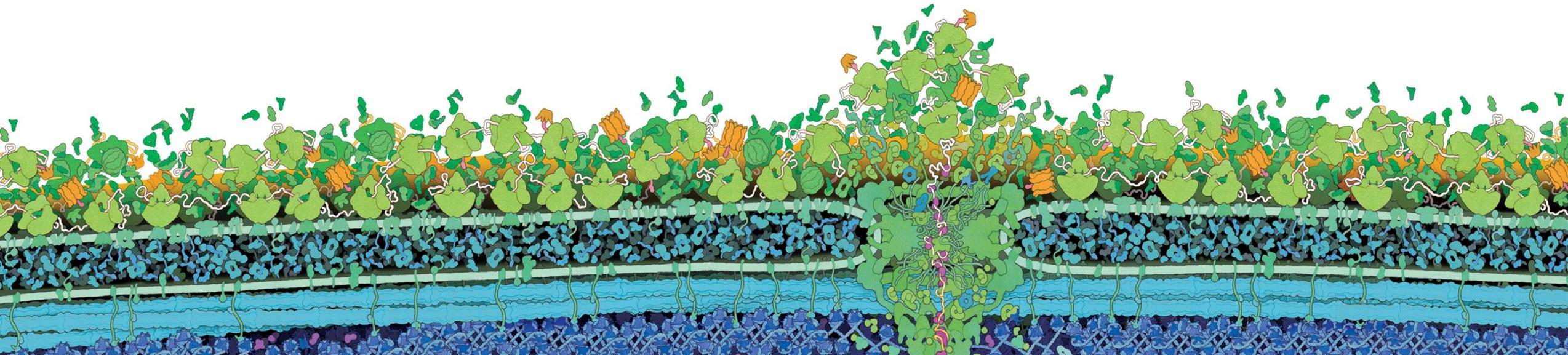
### Specificity



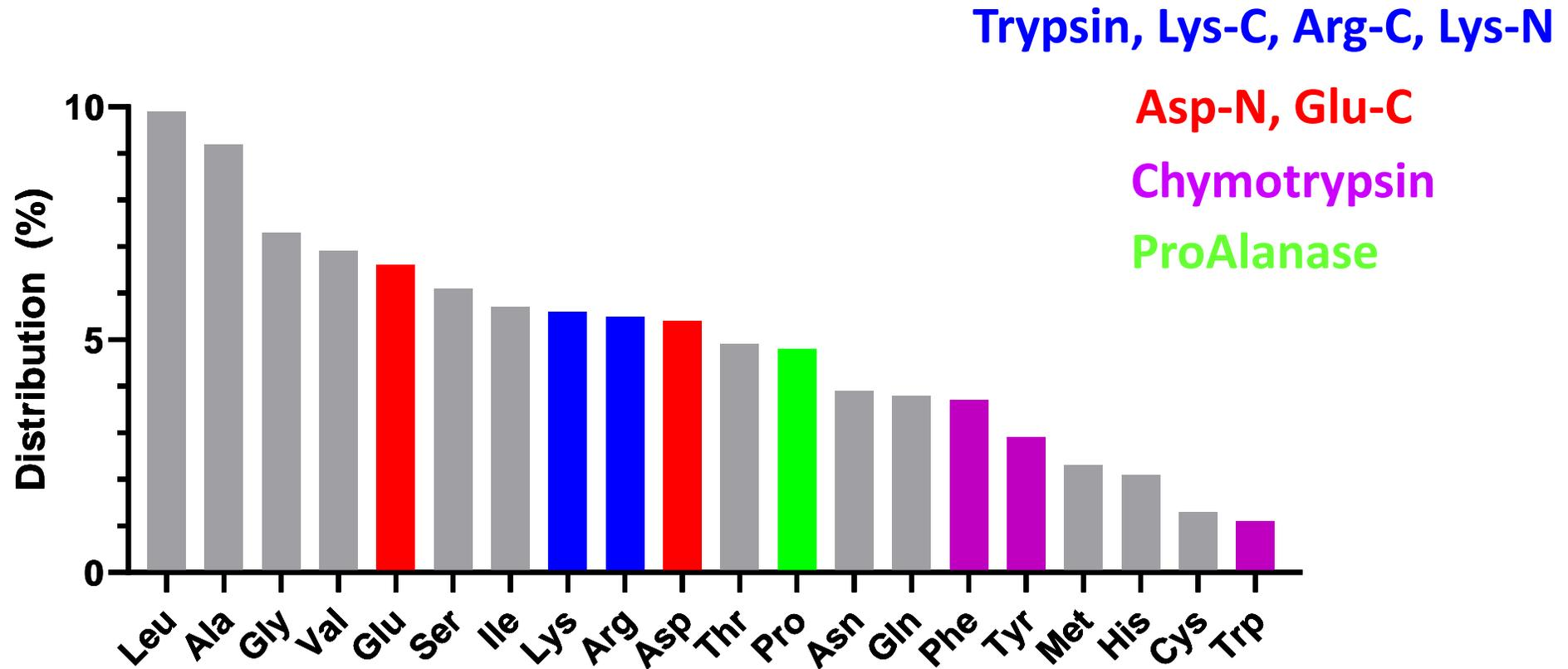
Cleavage efficiency at Aspartic acid improves slightly  
Increased # of cleavages at Glutamic acid reduces specificity

# ProAlanase

New Protease with Specificity for Proline and Alanine



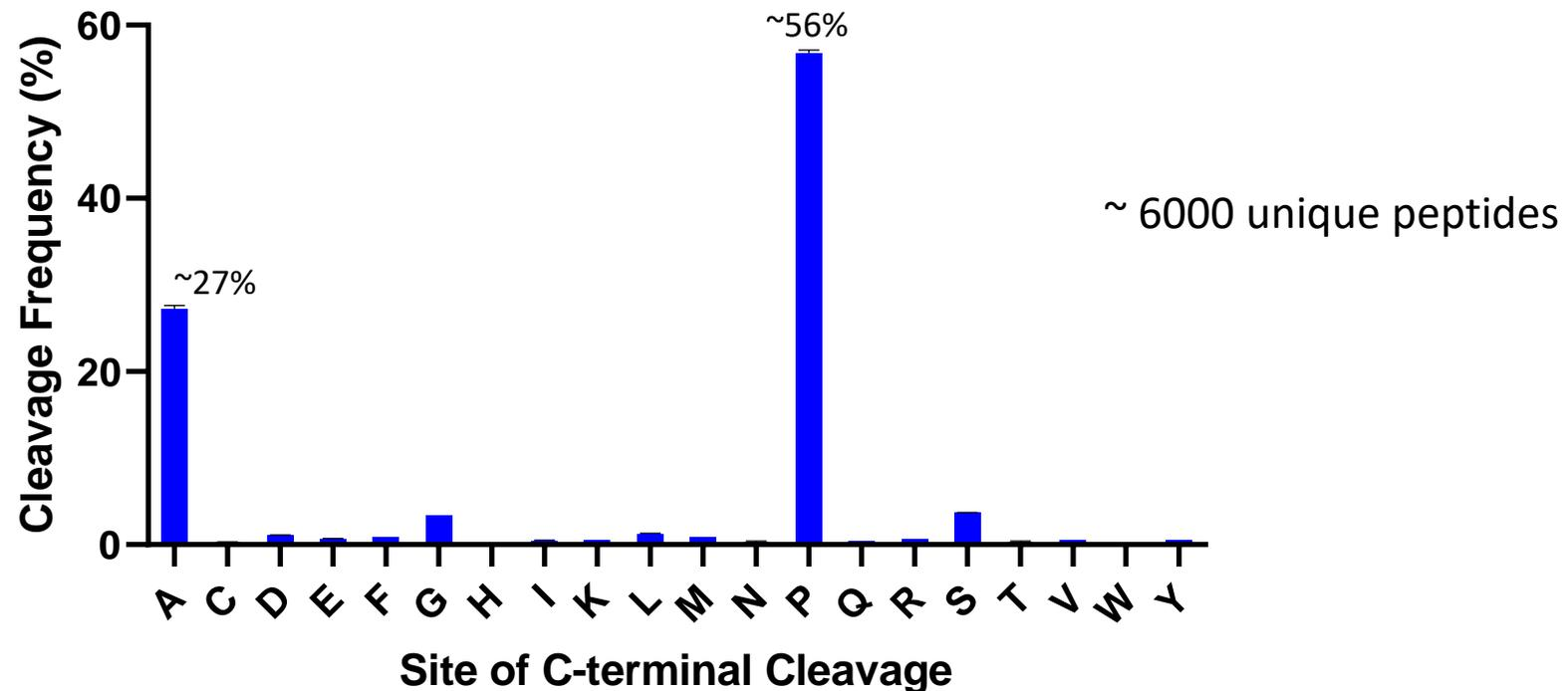
# Current Proteases do not Address the Full Proteome



Source: UniProt database, Sept 2019

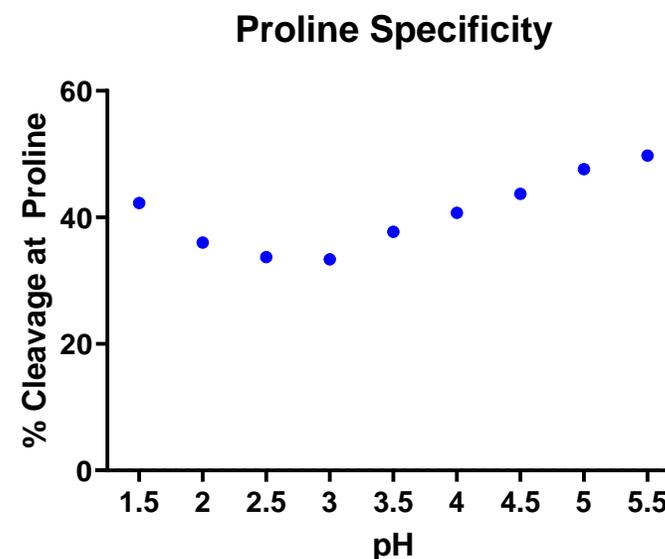
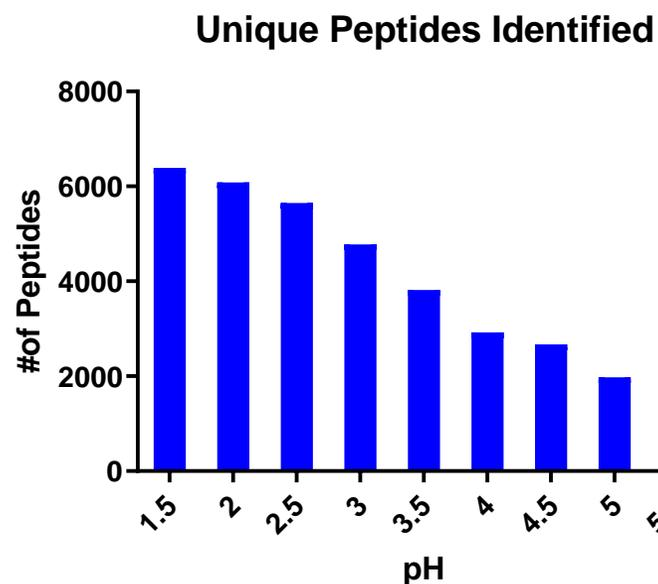
# ProAlanase Cleaves Primarily at Proline and Alanine

- ProAlanase is a serine protease from *Aspergillus niger* active at low pH
- Human K562 extract was digested with ProAlanase at pH 1.5 for 2 hours at 37°C at a 1:100 E:S ratio.
- Data collected with a Q Exactive Plus.
- Data were searched with Byonic (Protein Metrics) with no enzyme specified.



# ProAlanase Performance is Optimal around pH 1.5

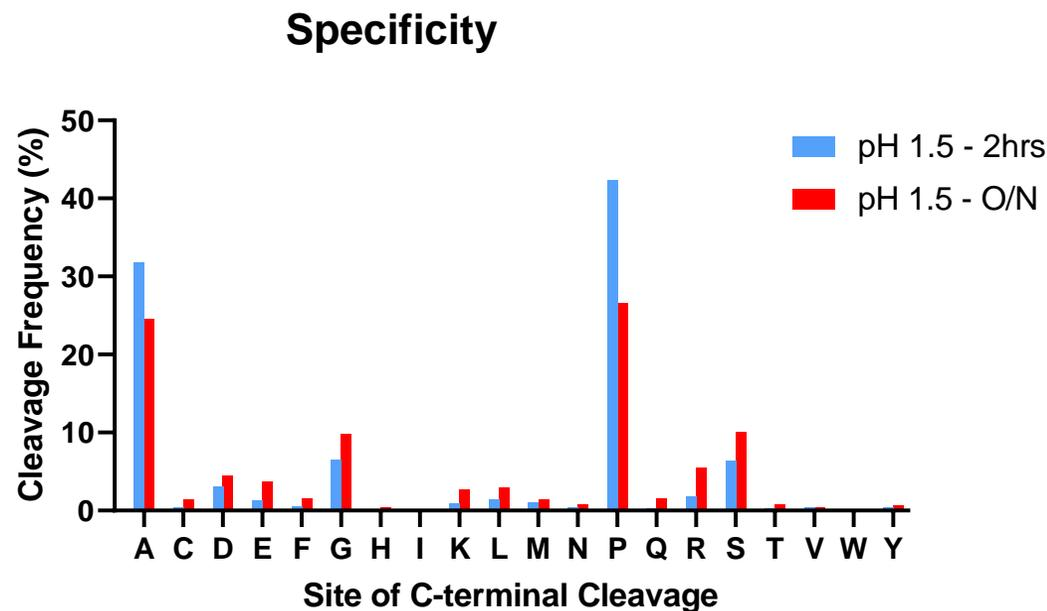
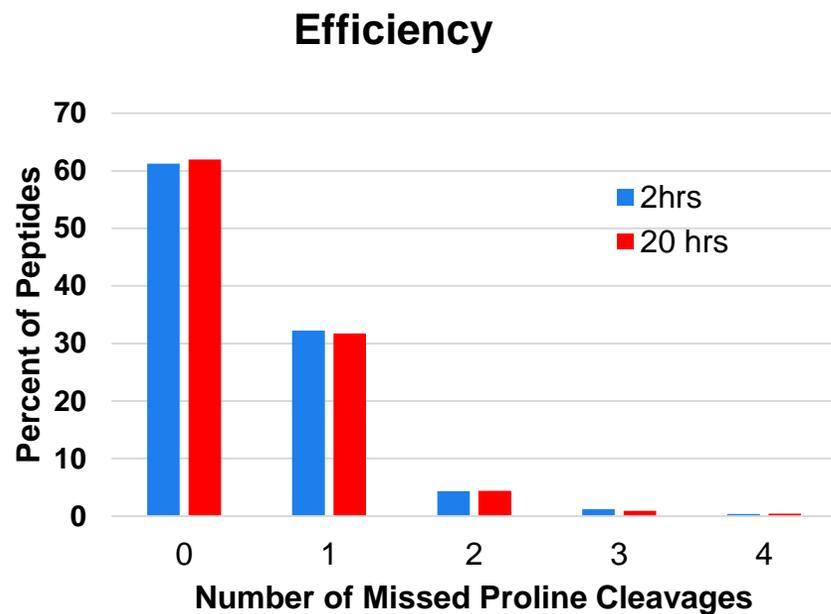
- Human K562 extract digested with ProAlanase *at various pH values* for 2 hours at 37°C at a 1:50 E:S ratio.



Best combination of specificity and peptide IDs observed at pH 1.5

# ProAlanase Performance is Optimal with Short Digestions

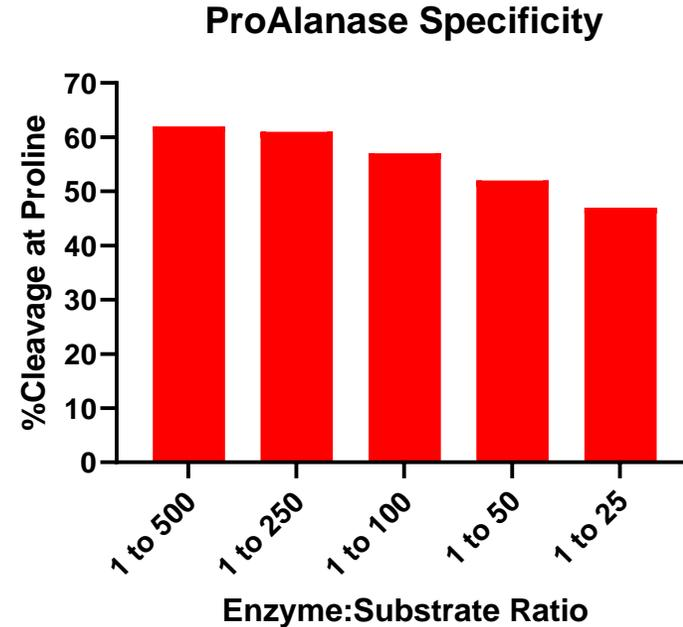
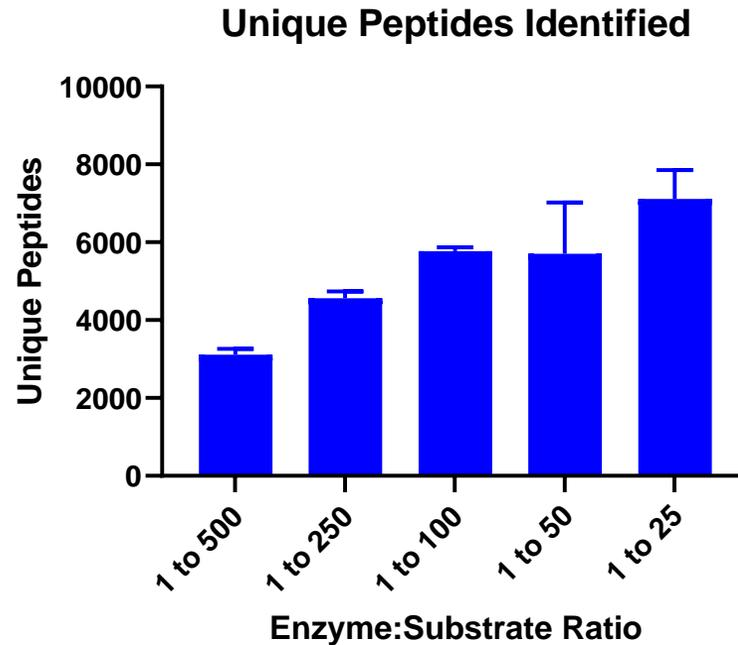
- Human K562 extract digested with ProAlanase at pH 1.5 *for 2 hours or overnight* at 37°C at a 1:50 E:S ratio.



Increasing digestion time beyond 2 hours does not increase efficiency but reduces specificity

# Changing ProAlanase Amount Affects Performance

- Human K562 extract digested with ProAlanase *at various E:S ratios* for 2 hours at 37°C at pH 1.5.



Increasing ProAlanase increases peptides IDs but reduces specificity

# How to Terminate Digestions with ProAlanase?

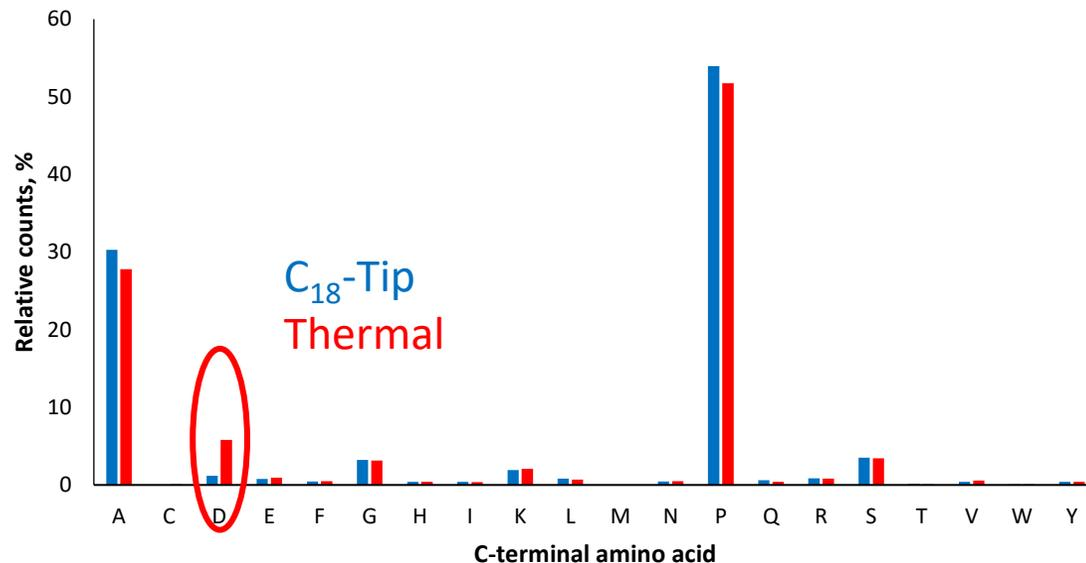


- Acidification with TFA or Formic Acid does not inhibit ProAlanase
  - Over-digestion of samples will occur unless reaction is terminated

Option 1: Heat 10 minutes at 90-95°C

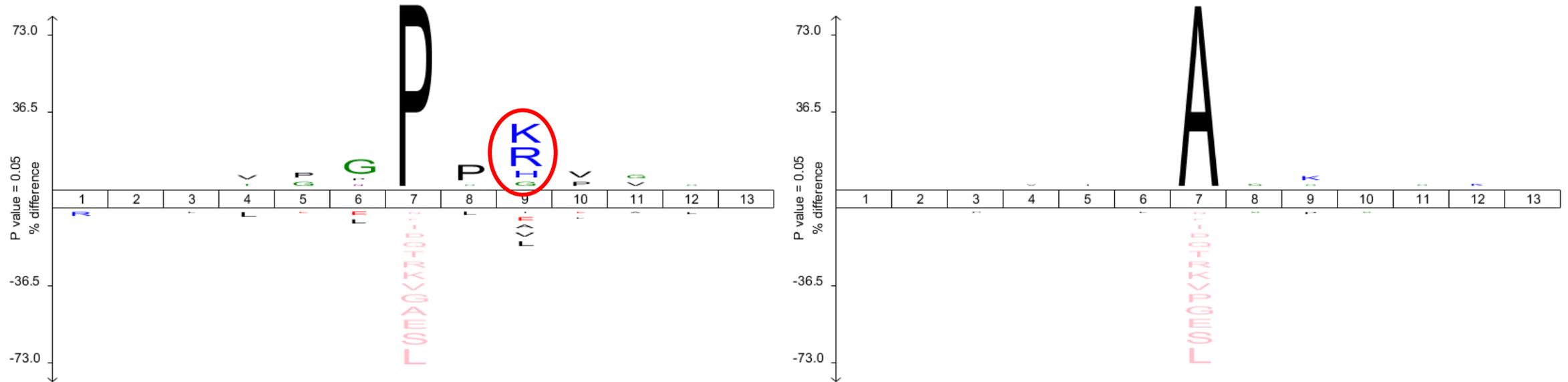
Option 2: C<sub>18</sub>-based cleanup (e.g. C<sub>18</sub> Tips)

Effect of Termination Method on Specificity



Termination Method	Ease of Use	Fragmentation?
Heat-based	Easy	Yes: D-P bonds
C <sub>18</sub> cleanup	Cumbersome (multiple samples)	No

# IceLogo Analysis of ProAlanase Cleavage Patterns

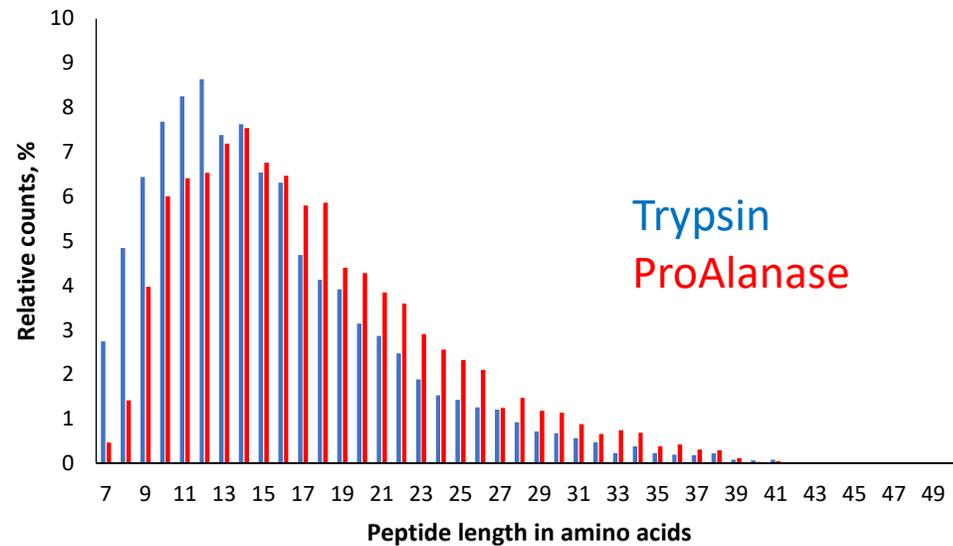


- Missed proline cleavage typically due to basic residue in P2' position.
- No clear patterns on missed cleavage at alanine

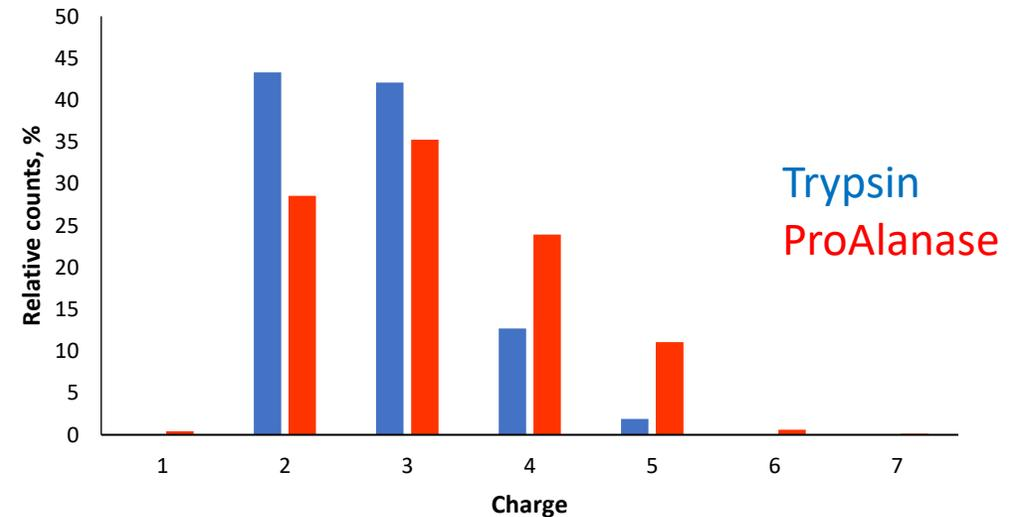
# Characteristics of ProAlanase-Derived Peptides



### Peptide Length Distribution

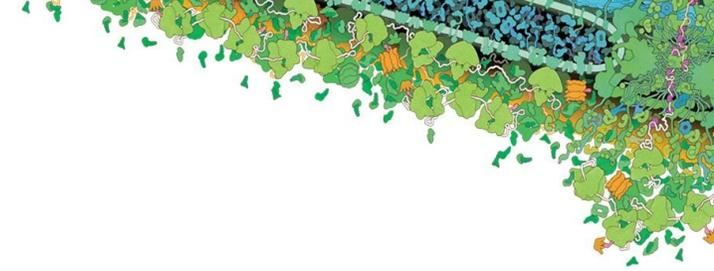


### Charge State Distribution

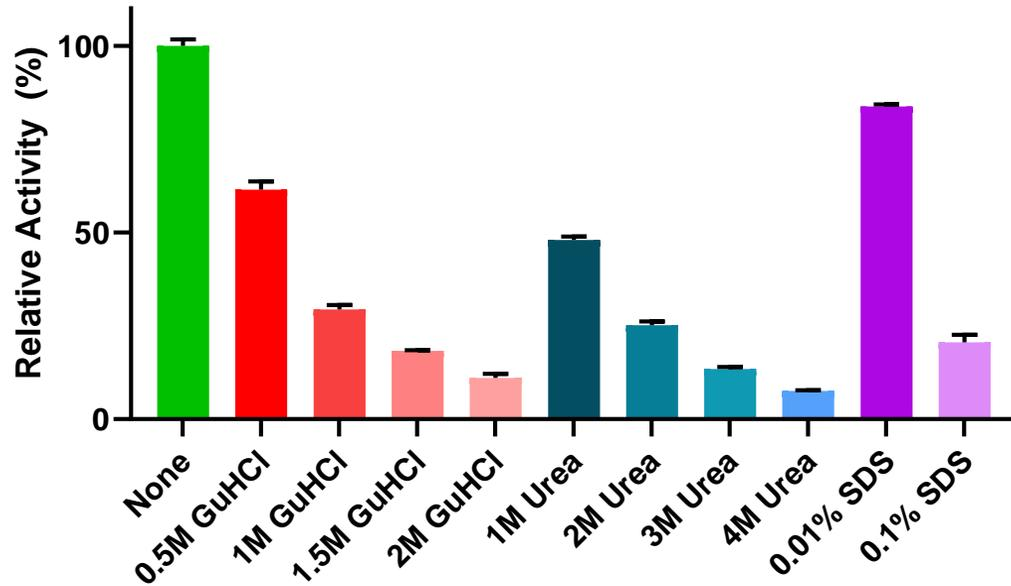


- Longer peptides than produced from trypsin
- More peptides at higher charge states than produced from trypsin

# Effect of Denaturants on ProAlanase Activity

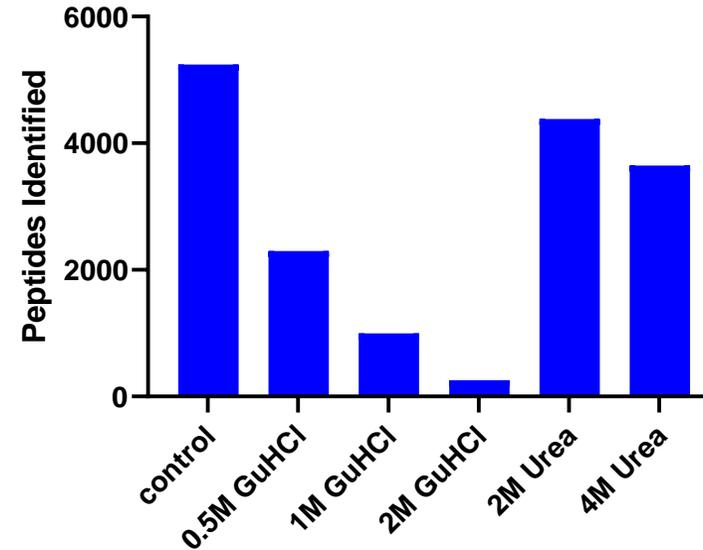


### Effect of Denaturants on ProAlanase Activity



### Synthetic Peptide Fluorescent Assay

### Effect of Denaturants on K562 Digestion with ProAlanase

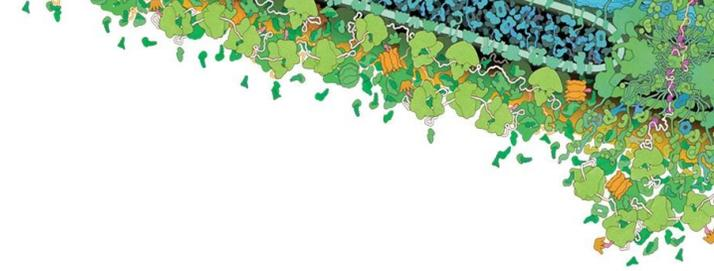


### K562 Digest & LC-MS/MS

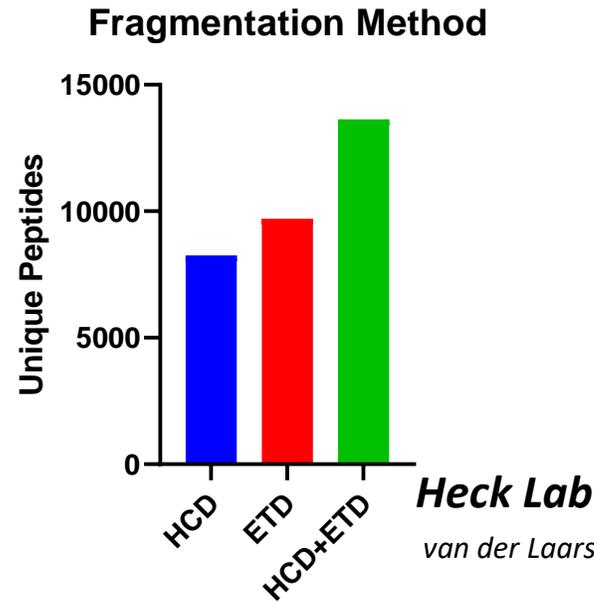
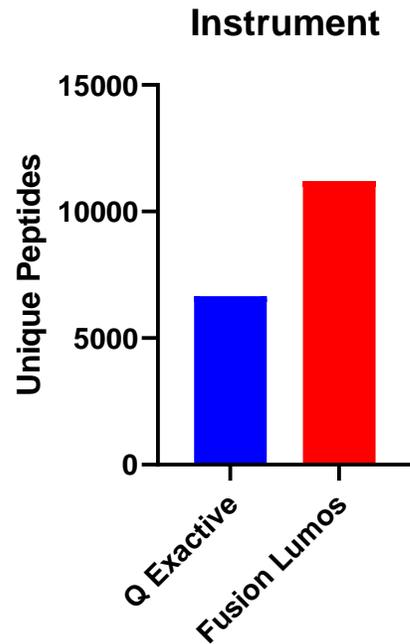


- 2-4M urea preferred if denaturant is needed
- \*Additional denaturants may not be needed at pH 1.5!

# MS Instrument and Software Affect Results

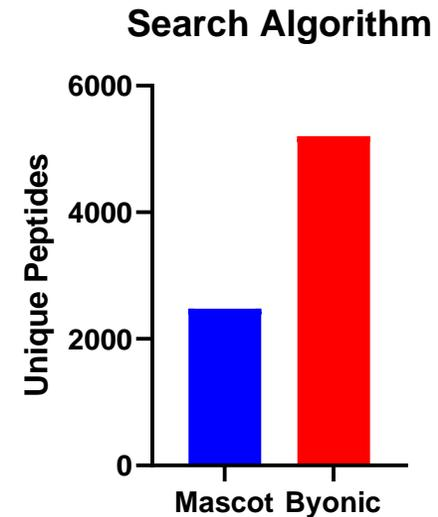


## Instrumentation



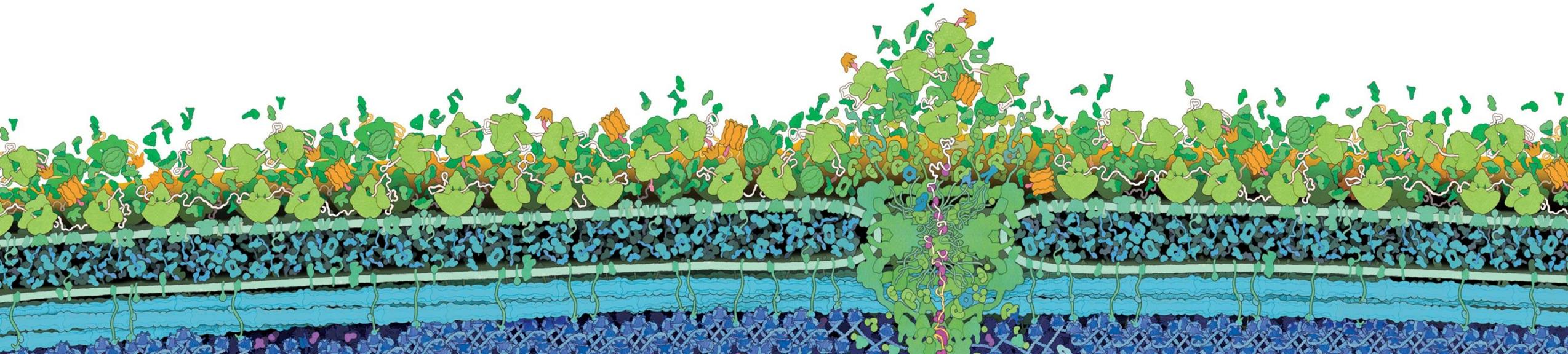
**Heck Lab**  
*van der Laarse et al (2019)*

## Software



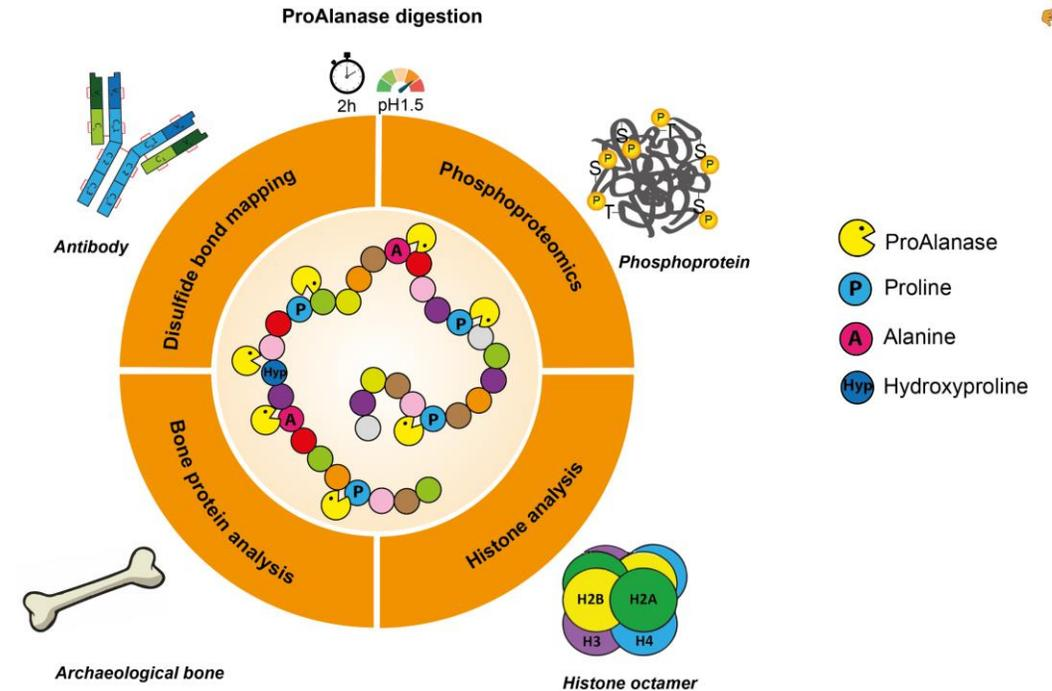
Optimize data collection and search strategies to increase peptide IDs

# Applications



# Applications of ProAlanase

- Proteomics
- Protein characterization / Peptide mapping
- Disulfide bond mapping
- *De novo* sequencing
- Histone characterization
- HDX-MS



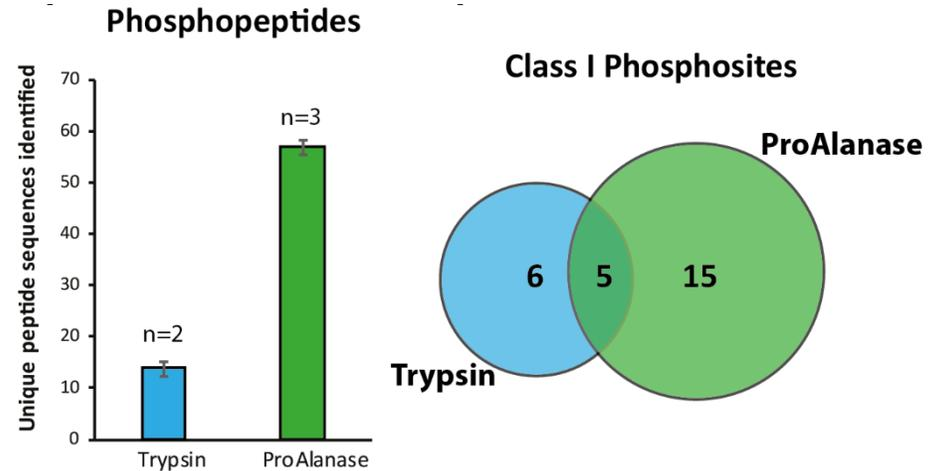
## New: Publication in Press

Samodova *et al.*, ProAlanase is an effective alternative to trypsin for proteomics applications and disulfide bond mapping.  
<https://www.mcponline.org/content/early/2020/10/05/mcp.TIR120.002129>

# Increase Sequence Coverage and Phosphosite Mapping



Data generated from in-gel digestion after I.P.

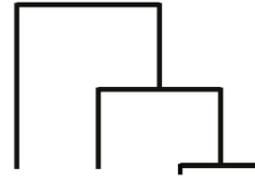


Increased sequence coverage  
 Increased phosphosite IDs

# Histone Characterization



## Pfam enrichment analysis

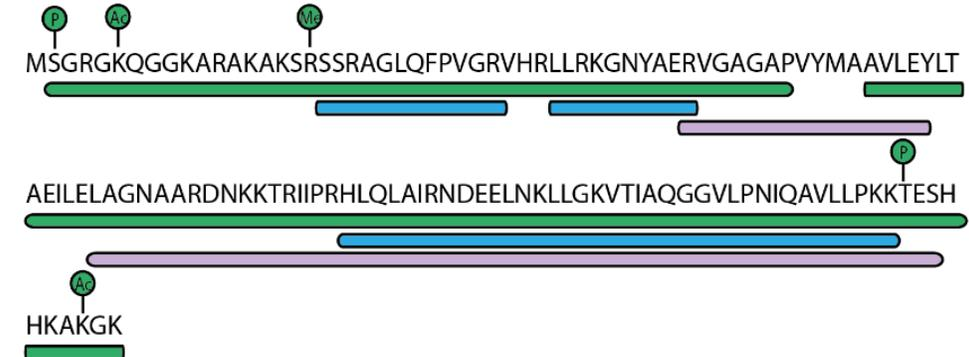


Term ID	Pfam protein domain (FDR-filtered)	Trypsin	Asp-N	Glu-C	ProAlanase
PF08534	Redoxin	Low	Low	Low	Low
PF10417	C-terminal domain of 1-Cys peroxiredoxin	Low	Low	Low	Low
PF16211	C-terminus of histone H2A	High	High	High	High
PF00578	AhpC/TSA family	Low	Low	Low	Low
PF01111	Cyclin-dependent kinase regulatory subunit	Low	Low	Low	Low
PF00125	Core histone H2A/H2B/H3/H4	High	High	High	High
PF01023	S-100/IcaBP type calcium binding domain	Low	Low	Low	Low
PF01849	NAC domain	Low	Low	Low	Low
PF00085	Thioredoxin	Low	Low	Low	Low
PF13207	AAA domain	Low	Low	Low	Low
PF00406	Adenylate kinase	Low	Low	Low	Low
PF00808	Histone-like transcription factor (CBF/NF-Y) and archaeal histone	Low	Low	Low	Low
PF13848	Thioredoxin-like domain	Low	Low	Low	Low
PF00631	GGL domain	Low	Low	Low	Low
PF00244	14-3-3 protein	Low	Low	Low	Low
PF00091	Tubulin/FtsZ family, GTPase domain	Low	Low	Low	Low
PF03953	Tubulin C-terminal domain	Low	Low	Low	Low
PF00241	Cofilin/tropomyosin-type actin-binding protein	Low	Low	Low	Low

>sp|Q6FI13|H2A2A\_HUMAN

- ▬ ProAlanase, 91 % coverage
- ▬ Trypsin, 46 % coverage
- ▬ Glu-C, 30 % coverage
- Phosphorylation
- Acetylation
- Methylation

## Histone H2A type 2-A sequence coverage



- ProAlanase significantly improves coverage of Histones compared with other proteases
- Allows for characterization of important Histone PTMs

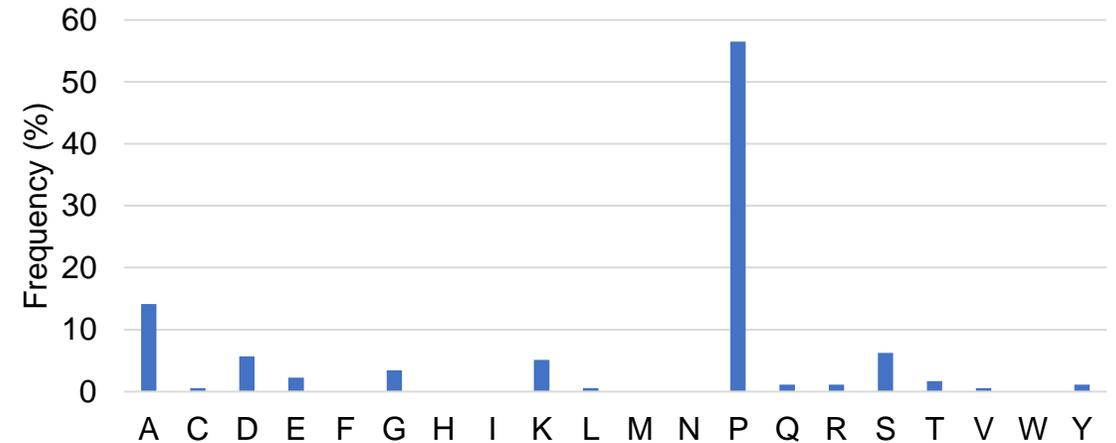
# Peptide Mapping of Biotherapeutic Proteins

- NISTmAb was digested with ProAlanaase at various E:S Ratios for 2 or 18 hours at 37°C

### Sequence Coverage

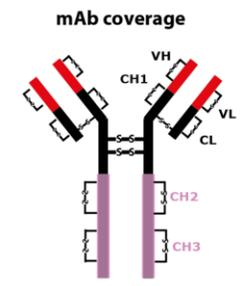
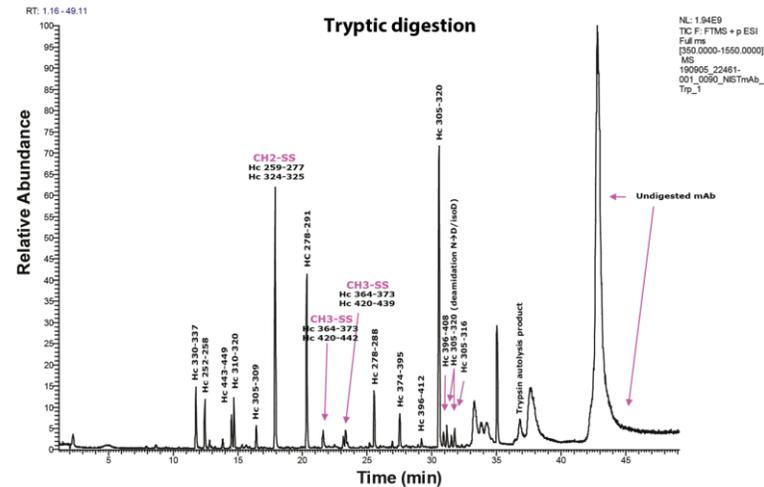
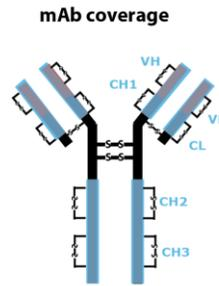
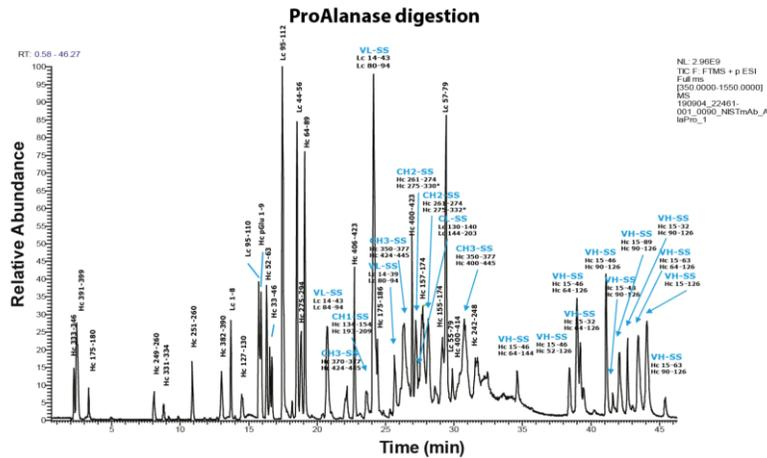
	1:250 E:S		1:100 E:S		1:50 E:S	
	1 Hr	18 Hrs	1 Hr	18 Hrs	1 Hr	18 Hrs
Heavy Chain	97	79	98	77	95	68
Light Chain	97	94	97	90	97	73

### Cleavage Specificity



- High coverage and high specificity obtained
- Short digestion times and low pH help minimize artifacts

# Disulfide Bond Mapping



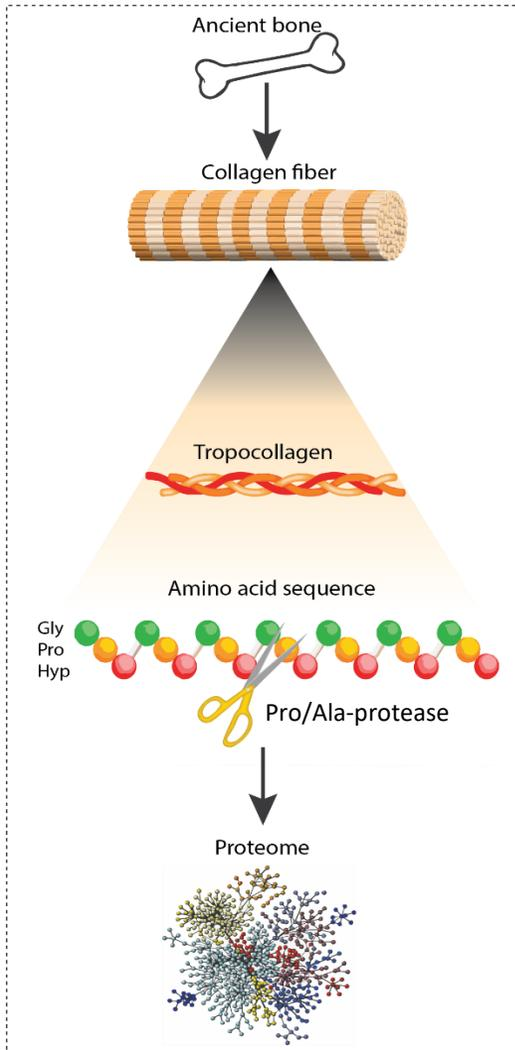
- Efficient digestion at pH 1.5
- Assignment of most S-S bonds

- Poor digestion at pH 8.0
- Assignment of few S-S bonds



- Low pH environment minimizes disulfide bond scrambling and artificial PTMs like deamidation
- Acidic pH acts as an effective denaturant of NISTmAb IgG allowing efficient digestion
- Assignment of all non-hinge disulfide bonds

# ProAlanase and Paleoproteomics



## Collagen

- Most abundant protein in bone (~ 90%)
- Well conserved thus not a good phylogenetic marker
- Suppresses other protein identification.
- Repetitive sequence motif: Gly-**Pro**-X & Gly-X-**Hyp**

**Aim:** Chop down collagen with ProAlanase and improve identification of other bone proteins.

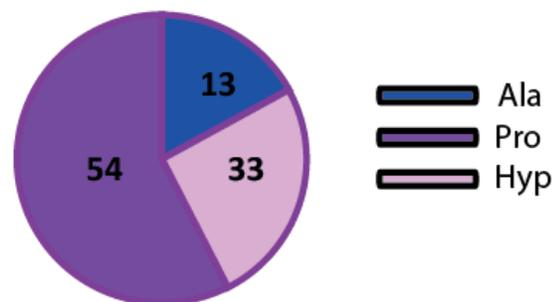
## Non-collagenous bone proteins can be used for:

- Reliable species identification
- Phylogenetic placement
- Disease identification

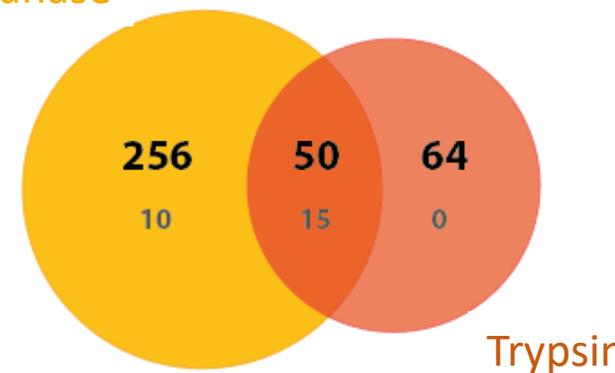
# Paleoproteomics of the Woolly Mammoth



ProAlanase Cleavage Sites in Bone



ProAlanase



*Protein groups IDd from mammoth bone*

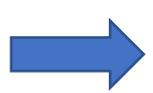
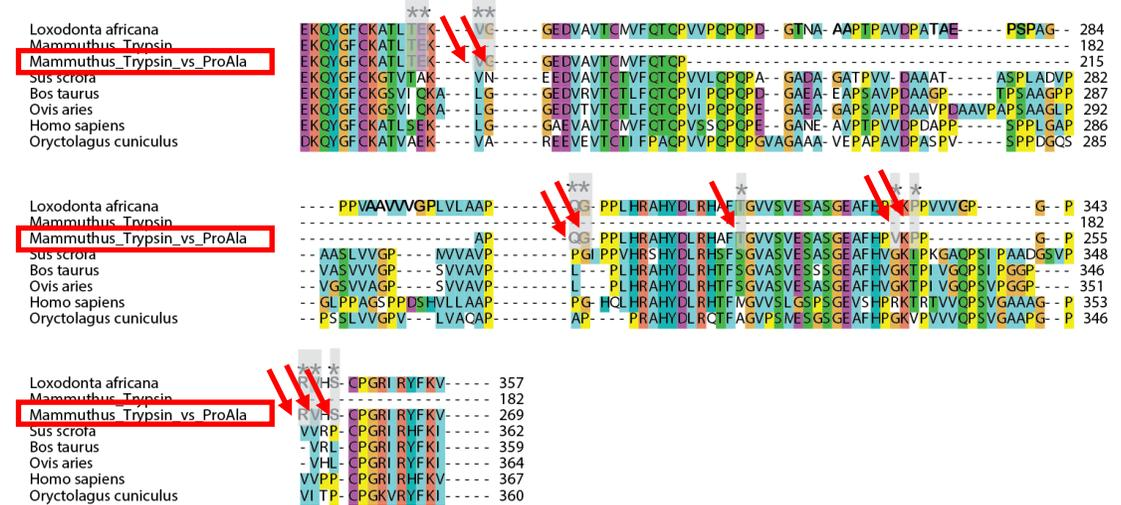
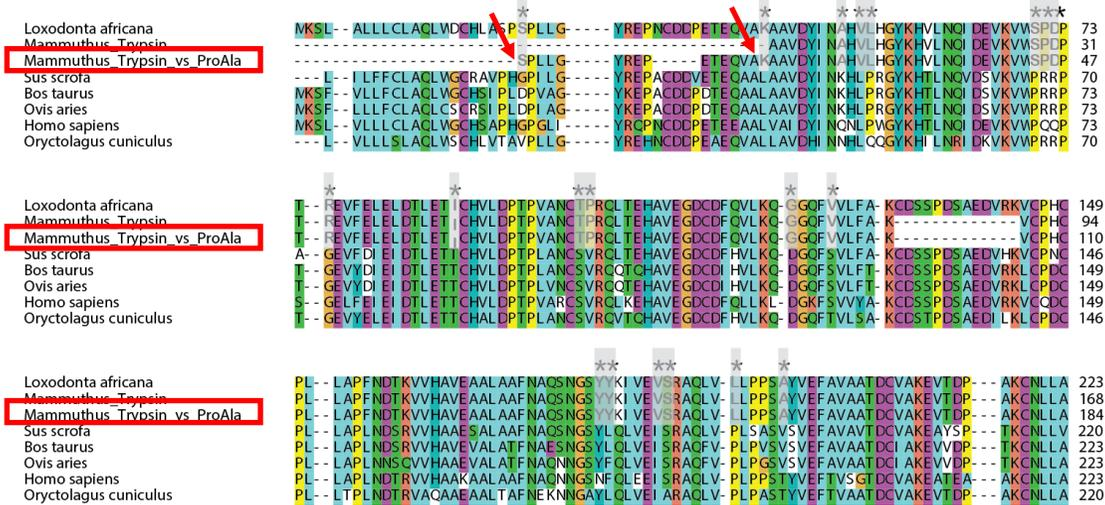


- ~3x more protein groups are identified in ProAlanase-digested sample
- 4x > complementary **non-collagenous** proteins (246), compared to trypsin (64)
- 10 complementary **collagen isoforms** identified with ProAlanase and 0 – with trypsin.

# Improved Sequence Coverage of Mammoth Fetuin



Fetuin-A sequence coverage across different species



- Combining peptides from trypsin and ProAlanase increases coverage of species-specific proteins.
- ProAlanase should be an effective tool for paleoproteomic analysis.

# ProAlanase Now Available

Product	Amount	Concentration
ProAlanase	5 µg	0.2 µg/µL
ProAlanase Plus	15 µg	0.5 µg/µL

## Ordering:

To purchase ProAlanase through our Early Access program, please contact your Promega sales representative directly for a quote. Alternatively, please email our early access team at [CAS@Promega.com](mailto:CAS@Promega.com)

## Technical Support:

For technical questions, please contact Chris Hosfield, Senior R&D Scientist [chris.hosfield@promega.com](mailto:chris.hosfield@promega.com)

# Acknowledgements

## University of Copenhagen

*Jesper Olsen*

*Diana Samodova*

Christian Kelstrup

Giulia Franciosa



Enrico Cappellini

## Novo Nordisk A/S

Christian Cramer

## Sapienza University of Rome

Maria Giuli

## Promega Corporation

*Mike Rosenblatt*

## Alpha Testers

Francis Crick Institute

Karolinska Institute

Lunenfeld-Tanenbaum Research Institute

Max Planck Institute of Biochemistry

ETH Zurich

University of Southern Denmark

EMD Serono

Amgen

MedImmune/AstraZeneca

MS Bioworks

Rapid Novor

Bioinformatics Solutions

Medical Proteoscope

JadeBio