

Examining the components of your peptide sample with AccuPep QC

Lauren Lu, Ph.D.

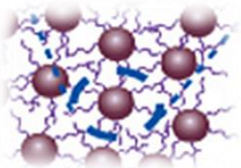
October 29, 2015, 9:00-10:00 AM EST



When do I need custom peptides?



Custom peptides play an important role in many research applications



Tissue Engineering

Hydrogels • stem cells
• wound healing



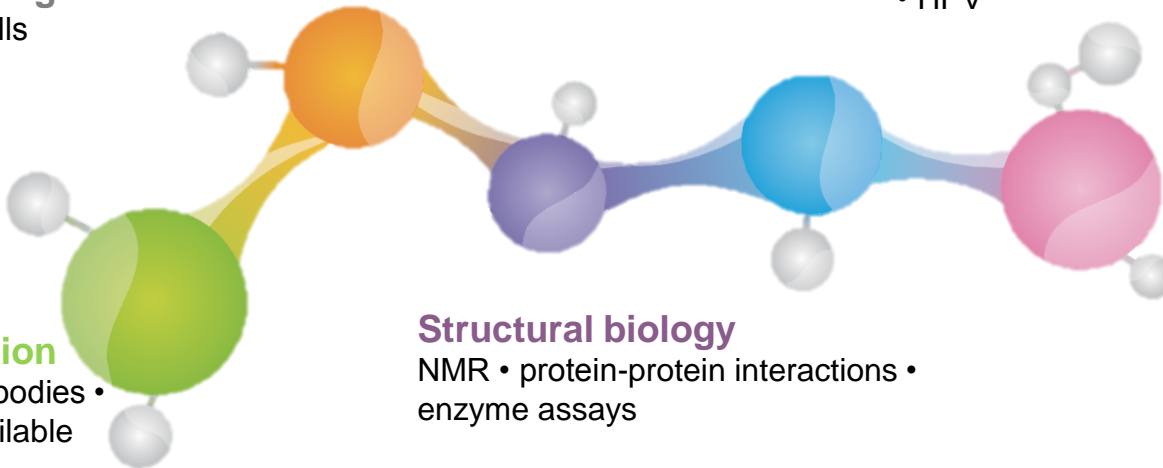
Drug Discovery

Antimicrobials • cancer • diabetes
• neurodegenerative diseases •
immunotherapy



Vaccine Development

HIV • cancer • influenza
• HPV



Antibody Generation

Phospho-specific antibodies •
non-commercially available
antibodies

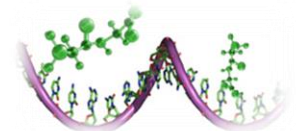


Structural biology

NMR • protein-protein interactions •
enzyme assays

Drug Discovery

siRNA



Common challenges with custom peptides



- ◆ Common issues that you might face
 - Unable to dissolve peptides with poor solubility
 - Improper dissolving way ruining peptide sample
 - Unknown contamination ruining your assay
 - Low experiment reproducibility from batch to batch
 - Non-reproducible results for quantitative experiments

How can these issues be addressed?



- ◆ For certain types of assays, additional testing is required to learn more about the contents of your peptide sample:
 - Appropriate solvents
 - Removal of TFA
 - The precise amount of net peptide
 - The presence of endotoxin
 - The water %
 - The pH value
 - The residual solvents
 - The identification of peptide impurities
 - ...

Benefits of analyzing your peptide content



GenScript offers a comprehensive QC service, AccuPep+, to help you get the most out of your custom peptides

- ◆ Reduces experimental troubleshooting
- ◆ Increases experiment reliability
- ◆ Ensure reproducible results

Features of the AccuPep+ Service



Quantification tests

- Amino Acid Analysis
- Peptide Content Analysis
- Counter-ion Quantification Analysis
- Moisture Content Analysis

Toxicity Tests

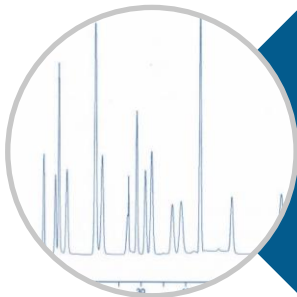
- TFA Removal and Analysis
- Endotoxin Analysis

Other tests

- Solubility Tests
- pH Test

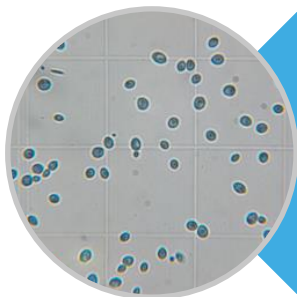
www.genscript.com/accupep_quality.html

AccuPep+ service test options



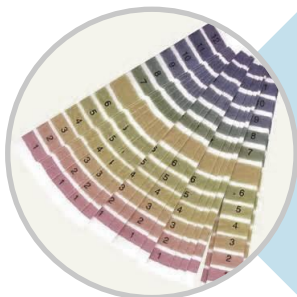
Quantification Tests:

- Do I really know all the possible components in my peptide sample?



Toxicity Tests:

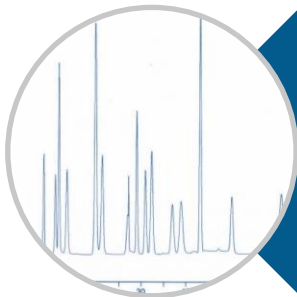
- What could make my experiment fail?



Other Tests:

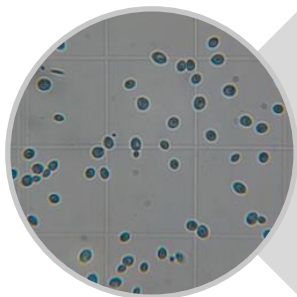
- What else can I do to accelerate my experiments?

AccuPep+ service test options



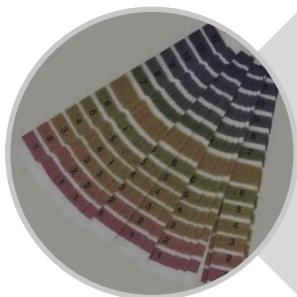
Quantification Tests:

- Do I really know all the possible components in my peptide sample?



Toxicity Tests:

- What could make my experiment fail?



Other Tests:

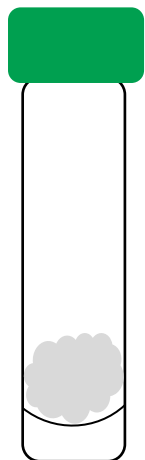
- What else can I do to accelerate my experiments?

Quantification Tests: Methods



Components

Quantification Methods



Target Peptide

Peptide Impurities

Amino Acid Analysis;
Peptide Content Analysis (Nitrogen
Element Analysis);
HPLC Analysis

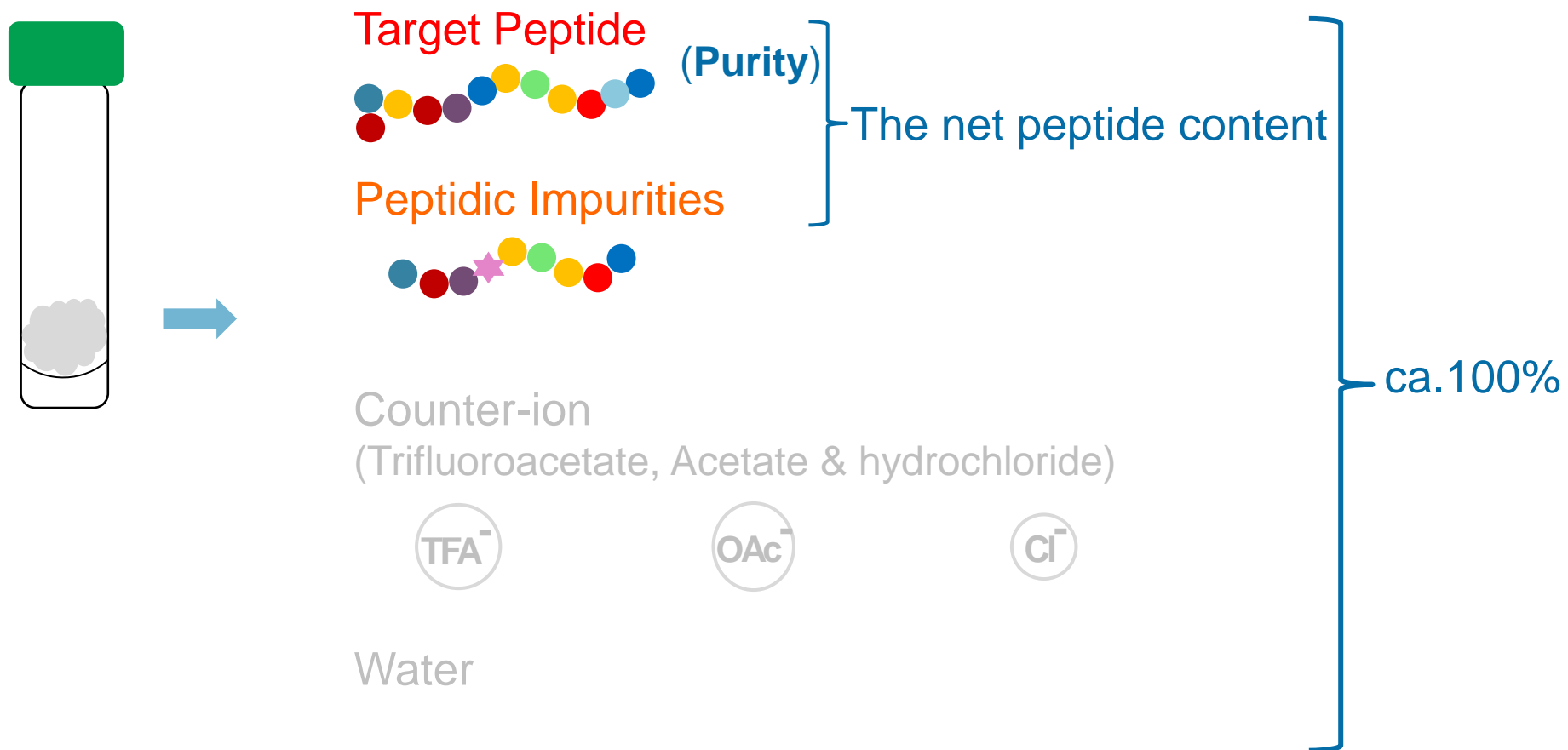
Counter-ion
(Trifluoroacetate, Acetate
& hydrochloride)

Counter Ion Quantification Analysis
(Ion chromatography)

Water

Moisture Content Analysis (Karl
Fischer coulometric titration)

How much peptide do I have?

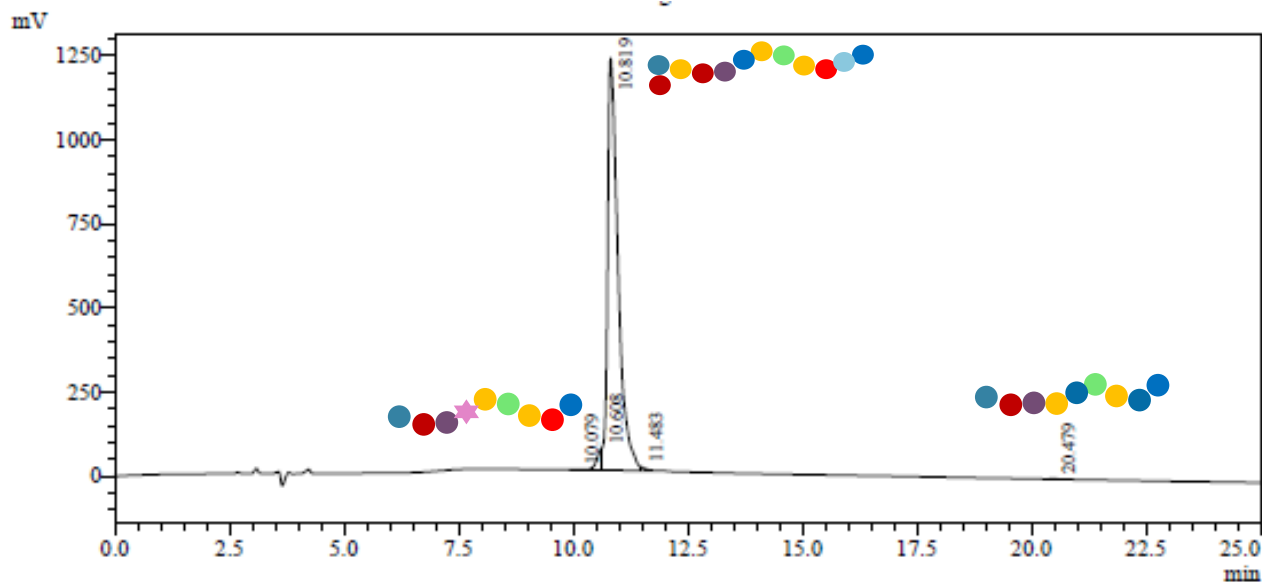


How is peptide purity measured?



◆ High performance liquid chromatography (HPLC) Purity

The ratio of peak area of target peptide in relation to all detected peak area



Detector A Ch1 220nm

Peak#	Ret. Time	Area	Height	Area %
1	10.079	12200	1255	0.061
2	10.608	441127	63265	2.201
3	10.819	19493777	1224629	97.257
4	11.483	83012	11085	0.414
5	20.479	13507	1763	0.067
Total		20043623	1301997	100.000

Peptide Purity: 97%

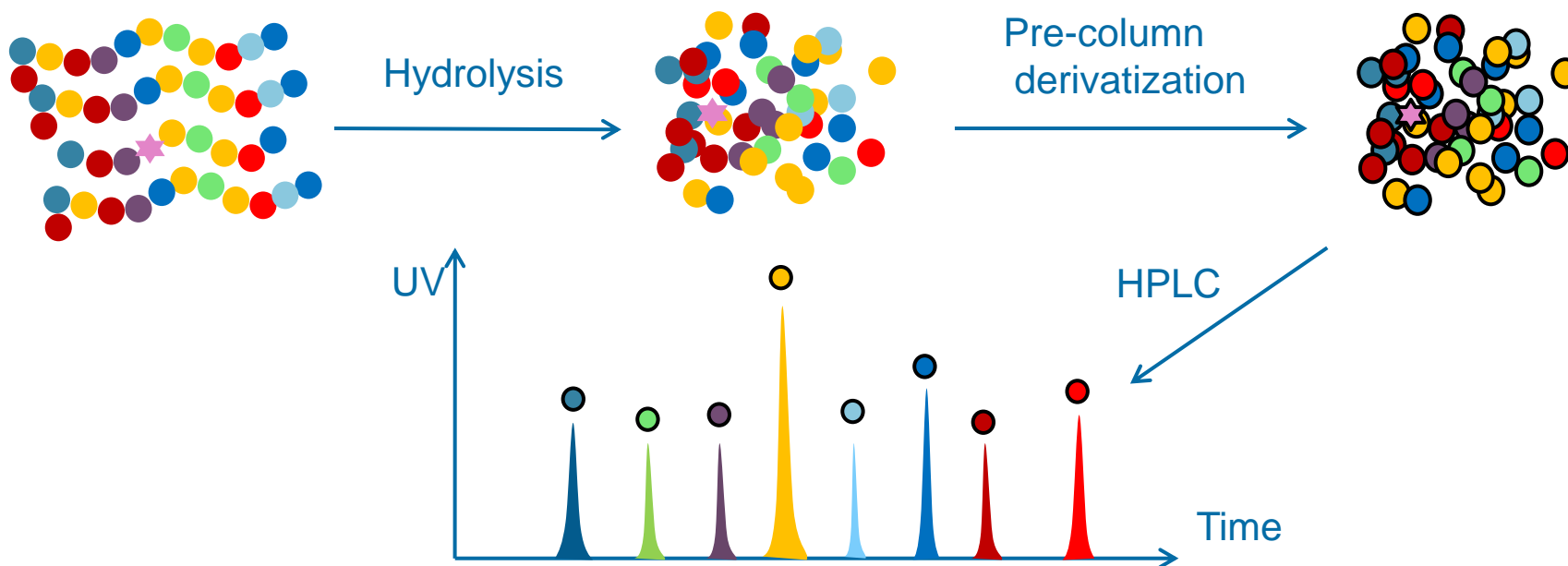
What is the amino acid composition of my peptide?



◆ Amino Acid Analysis (AAA) is ideal for:

- Determining the amino acid composition
- Determining the precise amount of net peptide in your sample

◆ Protocol:



Example AAA report



Peptide Sequence: VFNTRA

Amino acid (a.a.) residues	Theoretical a.a. number	The measured a.a. residues concentration ($\mu\text{mol/ml}$)	Measured a.a. number
Asp/Asn	1	0.1323	1.00
Glu/Gln			
Ser			
Gly			
His			
Arg	1	0.1308	1.00
Thr	1	0.1142	0.90
Ala	1	0.1245	1.00
Pro			
Tyr			
Val	1	0.1431	1.10
Met			
Cys			
Ile			
Leu			
Phe	1	0.1269	1.00
Trp			
Lys			

Tips:

1. Asn and Gln are de-aminated during hydrolysis to Asp and Glu.

2. Only highlighted a.a. residue is stable enough during hydrolysis to be used as for peptide content calculation.

Alternative way to determine the net peptide content



- ◆ **Peptide Content Analysis:** determines the precise amount of net peptide in the gross peptide sample
- ◆ Method
 - Nitrogen Element Analysis

Nitrogen %



$(C_x H_y N_z O_n S_m) \%$



Note: Given that the counter-ions (acetate, trifluoroacetate or hydrochloride) and the adsorbed water do not contain nitrogen, Nitrogen element analysis can be used for the net peptide content measurement.

How to calculate the net peptide



Delivered peptide

Delivered (Gross) Weight: 10 mg

Purity: 97%

Peptide content: 40%
(AAA)



Net peptide weight:

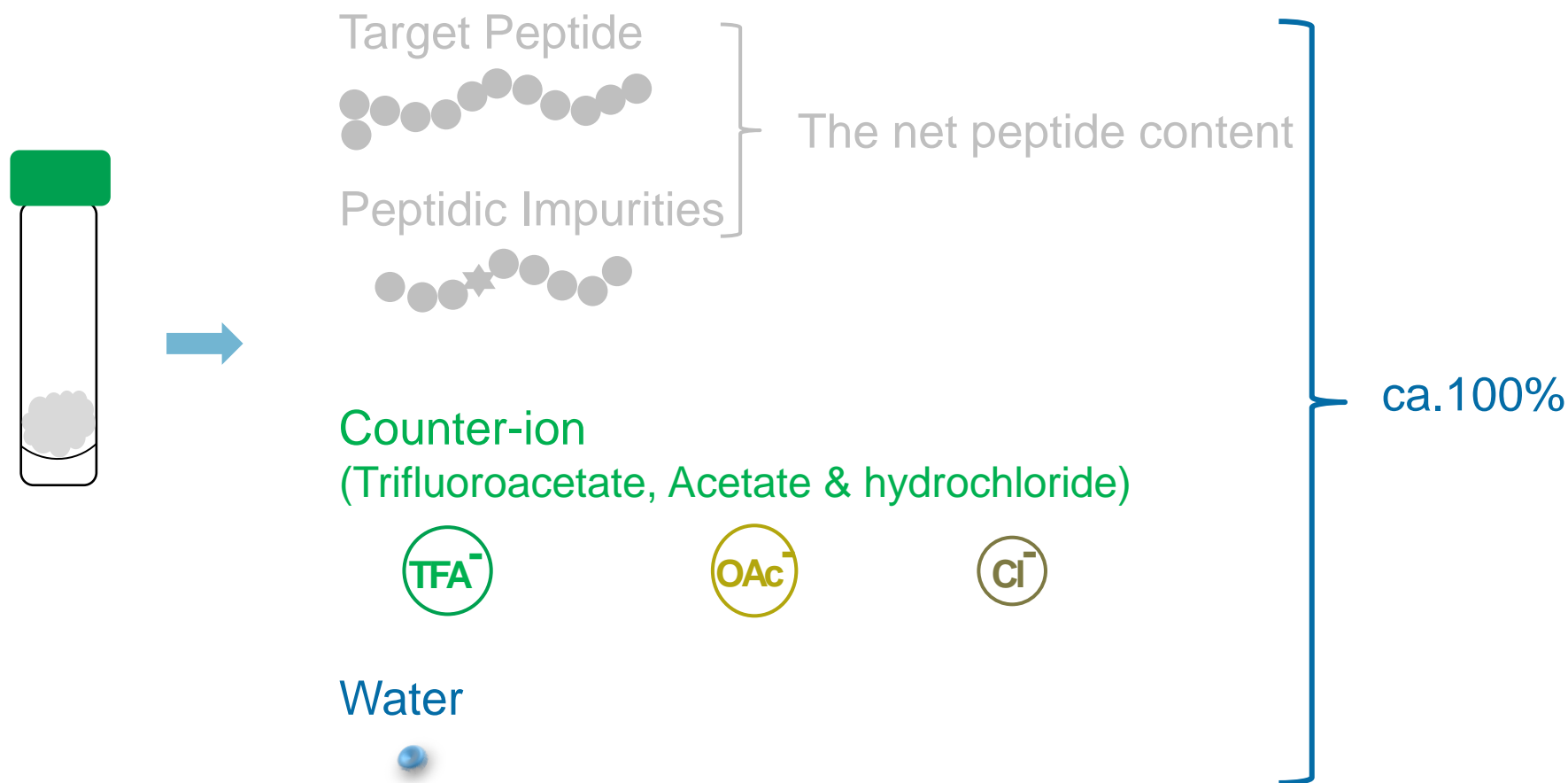
$$10 \text{ mg} * 40\% = 4.0 \text{ mg}$$

Exact amount of target peptide:

$$4.0 * 97\% = 3.88 \text{ mg}$$

With the purity and peptide content as determined by AAA, you can calculate the net peptide weight and exact amount of your peptide

Other constituents in your sample

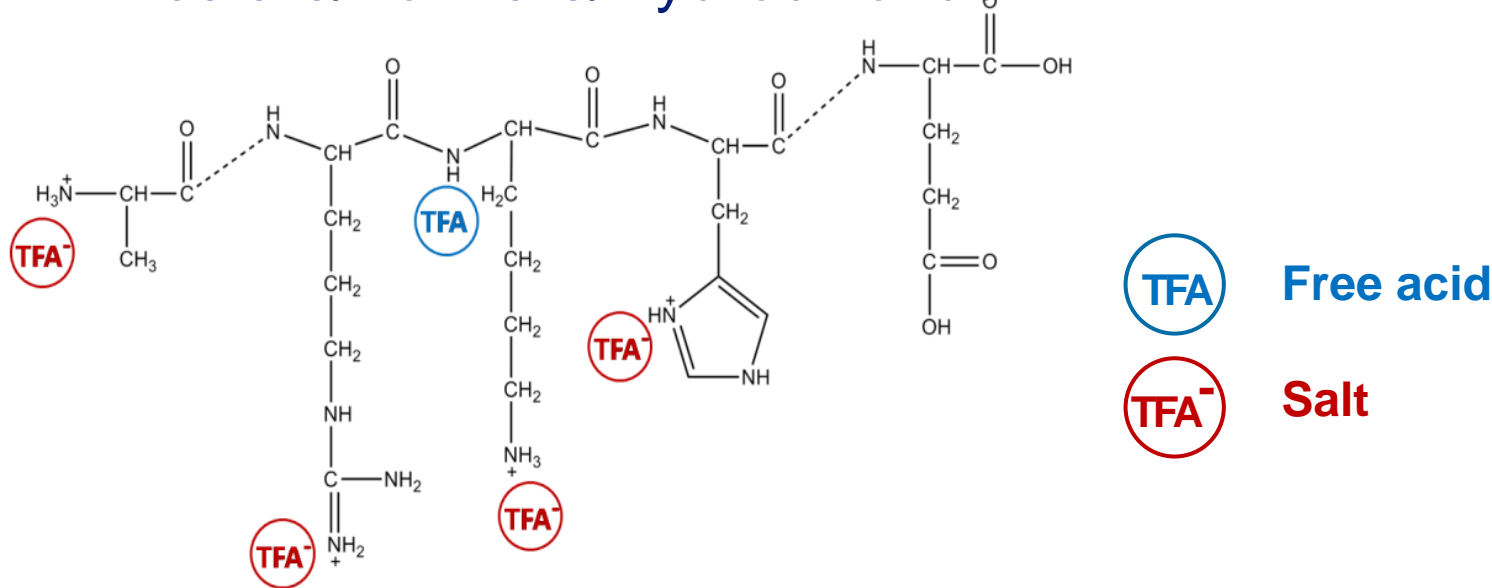


Counter-Ion Types and Sources



◆ Common counter-ion types

- **Trifluoroacetate** (**main type** for most delivered peptides)
- Acetate/Formate/Hydrochloride



◆ Counter-ion sources

- Peptide cleavage and purification
- Counter-ion exchange

Counter Ion Quantification Analysis

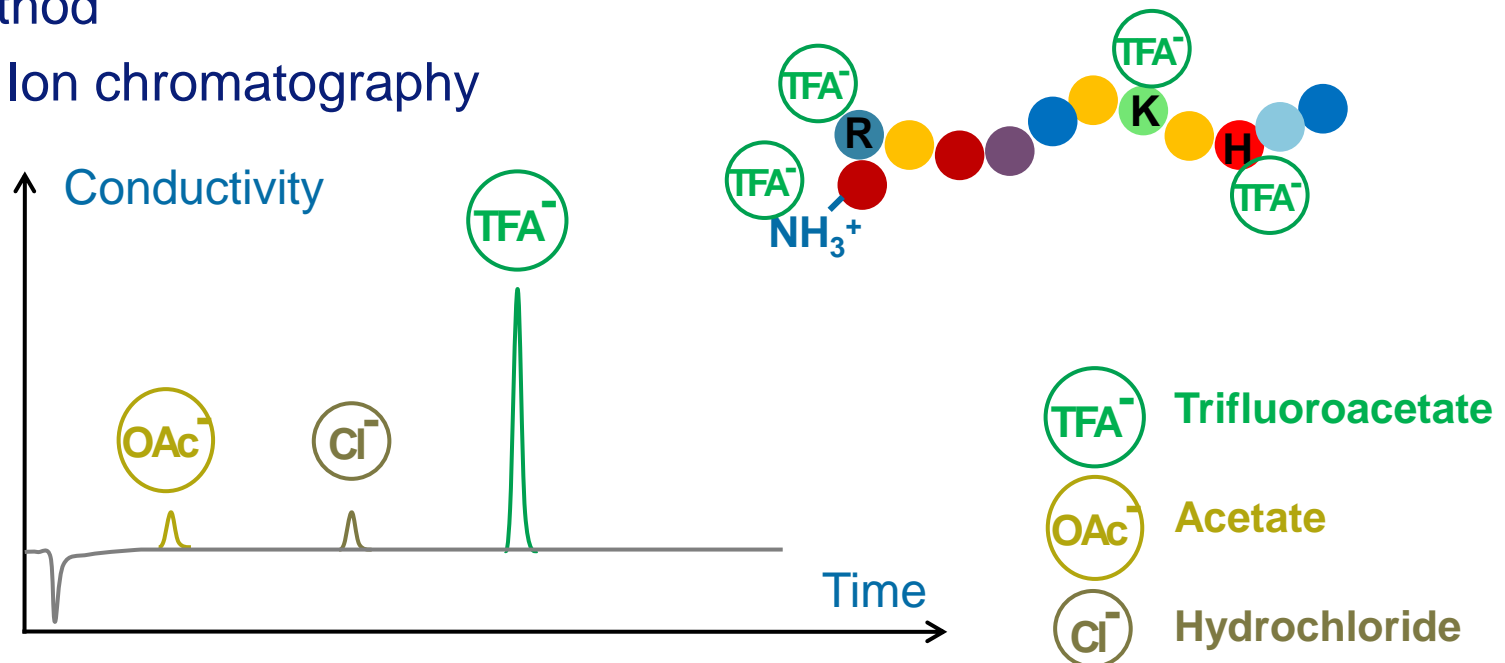


◆ Counter Ion Quantification Analysis is useful for:

- Cellular assays
- Active pharmaceutical ingredients (APIs)
- Manufactured products

◆ Method

- Ion chromatography



How can I estimate the theoretical TFA amount?



Peptide molecular weight: 1500 Da



TFA molecular weight: 114 Da



Peptide (adducted with counter-ion) molecular weight:
 $1500 + 114 \times 4 = 1956$ Da

The estimated TFA % = $114 \times 4 / 1956 = 23.3$ %

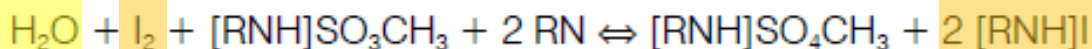
Moisture Content Analysis



- ◆ **Moisture Content Analysis** is useful for hydrophilic peptides that will retain the most water.

- ◆ **Method**

- Karl Fischer Coulometric Titration
- The basis:



I_2 reacts quantitatively with H_2O while the iodine is generated directly in the electrolyte by electrochemical means.

The rigorously quantitative relationship between the electric charge and the amount of iodine generated is used for high-precision dispensing of the iodine.





Components

Quantification Methods



Target Peptide

Peptide Impurities

Usually, 50-80%

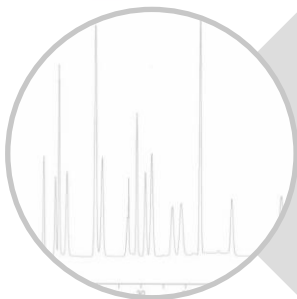
Counter-ion
(Trifluoroacetate, Acetate
& hydrochloride)

Usually, 30-10%
Dependent on basic a.a. residue
number

Water

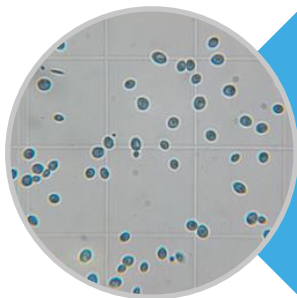
Usually < 10%,
but could be particularly high for
hydrophilic peptide

AccuPep+ service test options



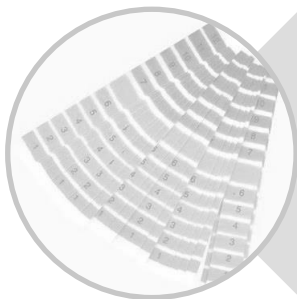
Quantification Tests:

- Do I really know all the possible components in my peptide sample?



Toxicity Tests:

- What could make my experiment fail?



Other Tests:

- What else can I do to accelerate my experiments?



◆ TFA removal and analysis

- TFA is a counter anion for normal peptides
- Trace amount of TFA can cause cytotoxicity in cell culture assays

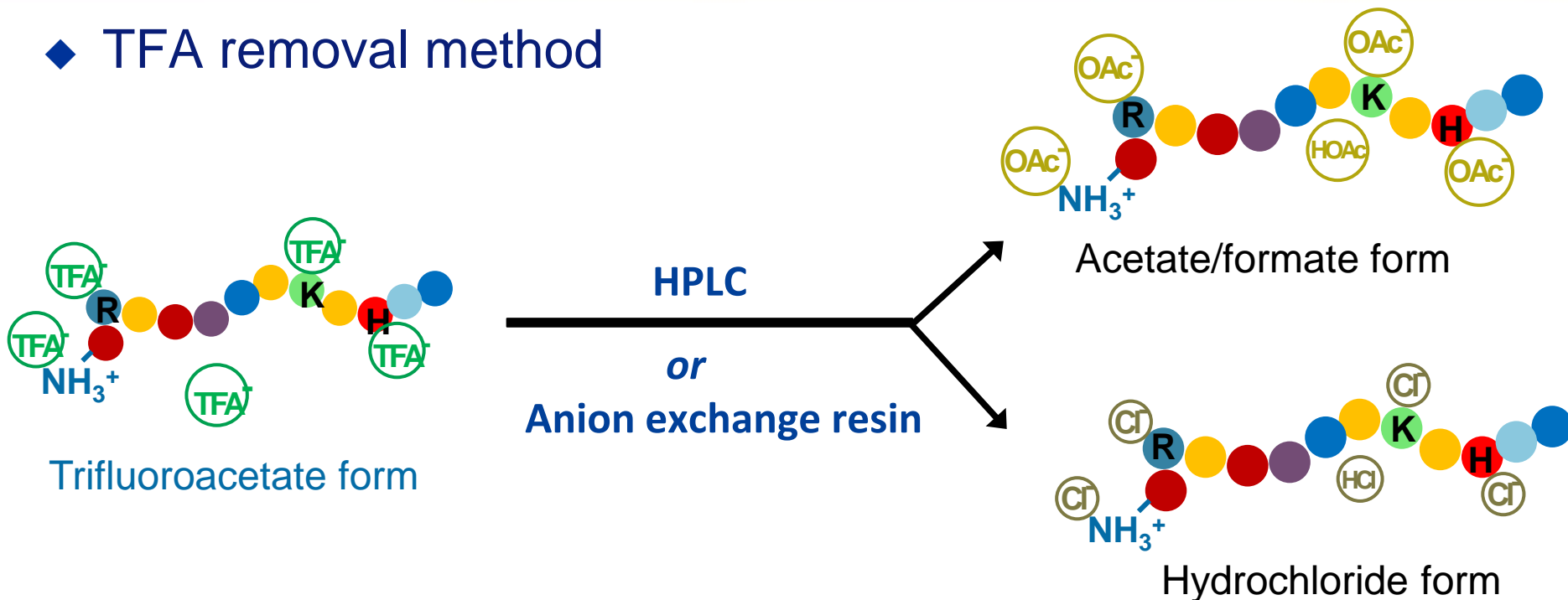
◆ Endotoxin Analysis

- Endotoxins are easily introduced into peptides during any process of peptide production.
- Small concentrations of endotoxin can decrease cell viability or cause immune response in cellular assays.

TFA Removal and Analysis



◆ TFA removal method



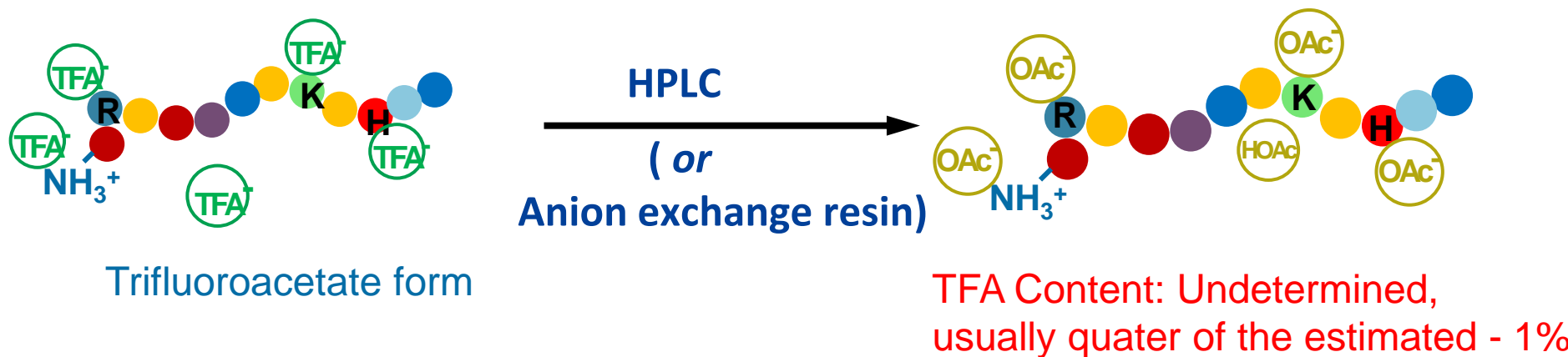
◆ How to choose counter-ion form (in the point view of production)

- OAc^- Suitable for unstable amino acids, such as Cys, Met, Gln at N-terminus
- Cl^- Suitable for peptides with low solubility

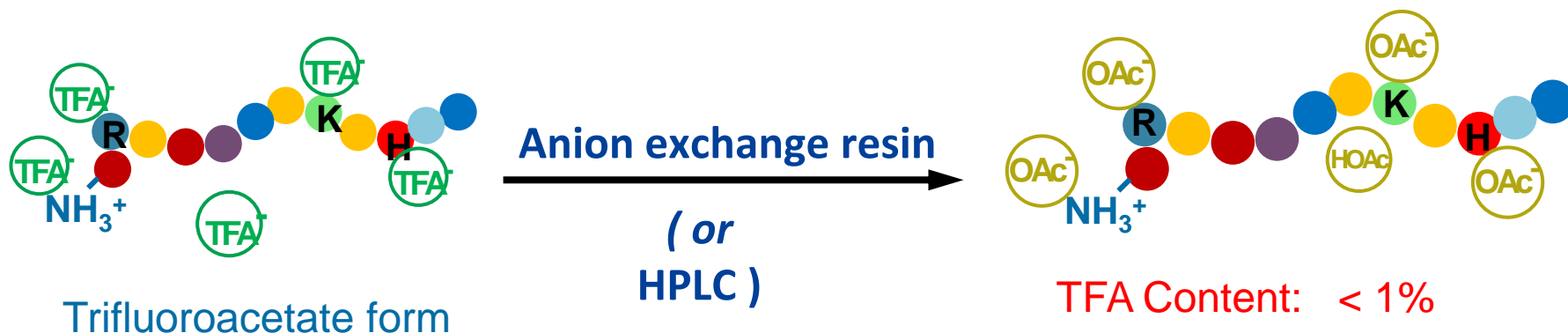
Two Types of TFA Removal Services



◆ Standard TFA Removal Service



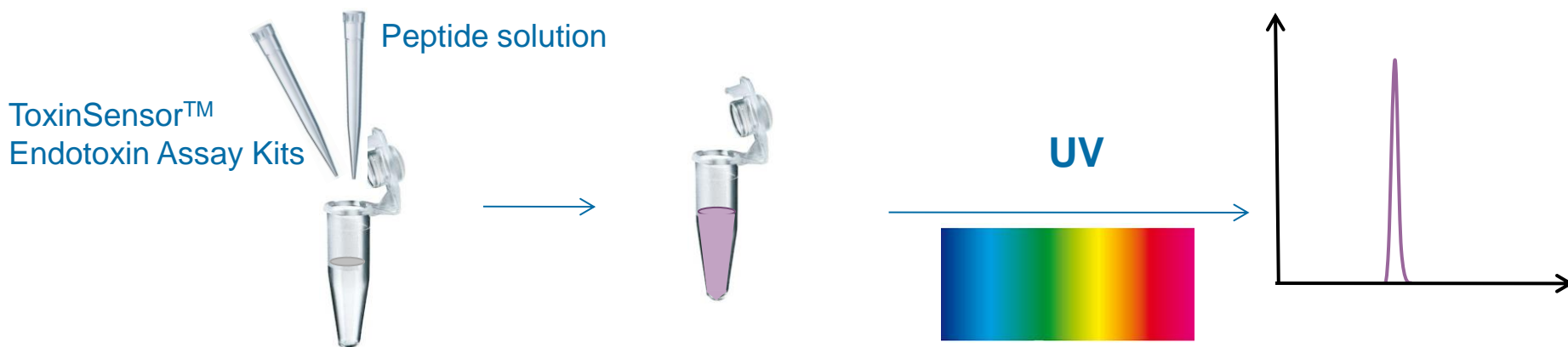
◆ Guaranteed TFA Removal Service



Endotoxin Analysis



- ◆ Endotoxins (lipopolysaccharides)
 - Major components of the cell walls of gram-negative bacteria
 - Introduced into custom peptides during peptide production
- ◆ Method
 - Chromogenic *Tachypleus* amebocyte lysate or *Limulus* amebocyte lysate test
 - Guaranteed high-sensitivity: 0.005 EU/ml



Which service is best for me?



- ◆ A variety of applications benefit from TFA removal or endotoxin analysis:

Application	Method
Analytical analyses that are influenced by TFA ions: <ul style="list-style-type: none">- Infrared (IR) spectroscopy- Circular dichroism (CD) spectroscopy	TFA removal and analysis
Cell culture assays	TFA removal and analysis
Cosmetics and pharmaceutical applications	TFA removal and analysis
Cell culture assays sensitive to endotoxin or prone to immune responses	Endotoxin analysis

Case study: effect of TFA on cell culture



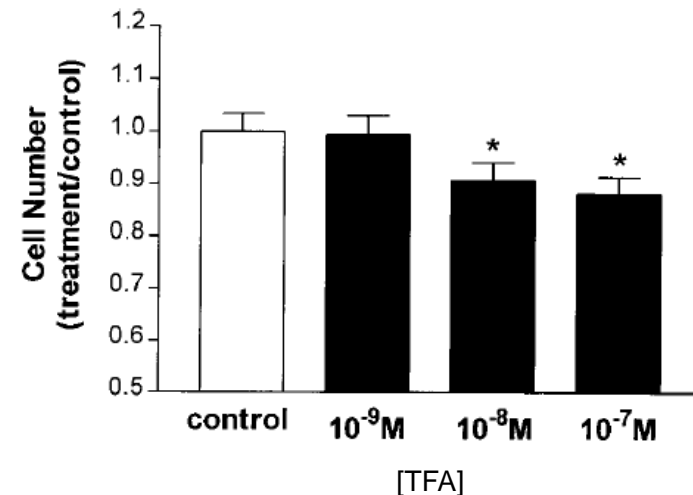
- ◆ Even trace amounts of TFA can cause cytotoxicity in cell culture assays

Trifluoroacetate, a contaminant in purified proteins, inhibits proliferation of osteoblasts and chondrocytes

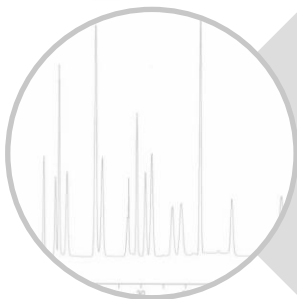
J. Cornish, K. E. Callon, C. Q.-X. Lin, C. L. Xiao, T. B. Mulvey, G. J. S. Cooper, I. R. Reid

American Journal of Physiology - Endocrinology and Metabolism Published 1 November 1999 Vol. 277 no. 5, E779-E783 DOI:

- ◆ Peptides containing TFA at concentrations ranging from 10^{-9} to 10^{-7} M were supplemented to osteocyte and bone cultures.
- ◆ Viability was assessed by [3 H]thymidine incorporation after 24 hours

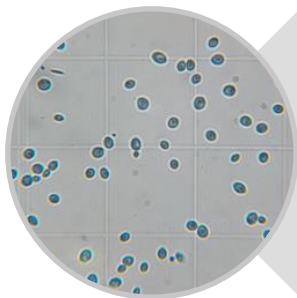


AccuPep+ service test options



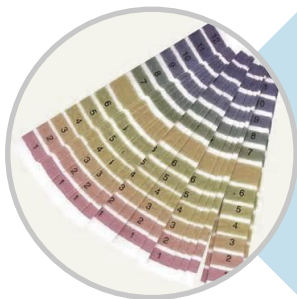
Quantification Tests:

- Do I really know all the possible components in my peptide sample?



Toxicity Tests:

- What could make my experiment fail?



Other Tests:

- What else can I do to accelerate my experiments?

Solubility Test

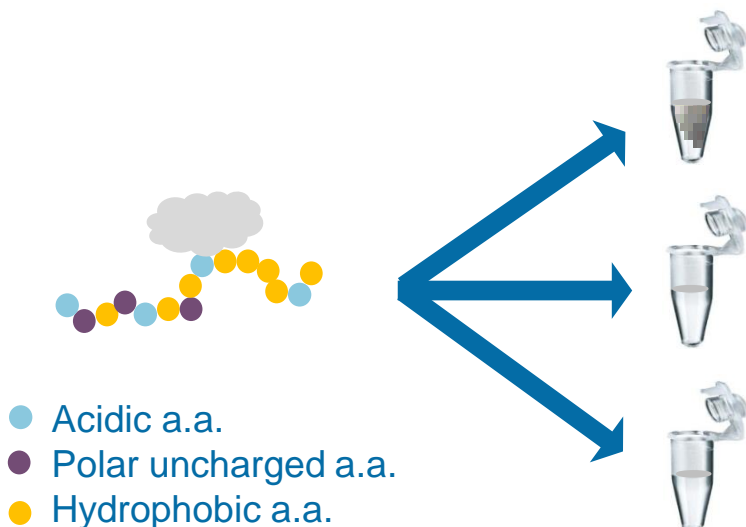


- ◆ Hydrophobic peptides
 - Containing > 50% hydrophobic amino acid
 - Length > 5 a.a.
- Hydrophobic :
 - Ala (A), Trp (W), Leu (L), Ile (I), Phe (F), Met (M), Val (V), Pro (P)
- Basic:
 - Arg (R), His (H), Lys (K)
- Polar uncharged:
 - Asn (N), Cys (C), Gly (G), Gln (Q), Ser (S), Thr (T), Tyr (Y)
- Acidic:
 - Asp (D), Glu (E)

Components of the solubility report



◆ What is included in your solubility report?



Benefits:

Reduces troubleshooting
Saves time and peptide products
Particularly useful for peptide libraries

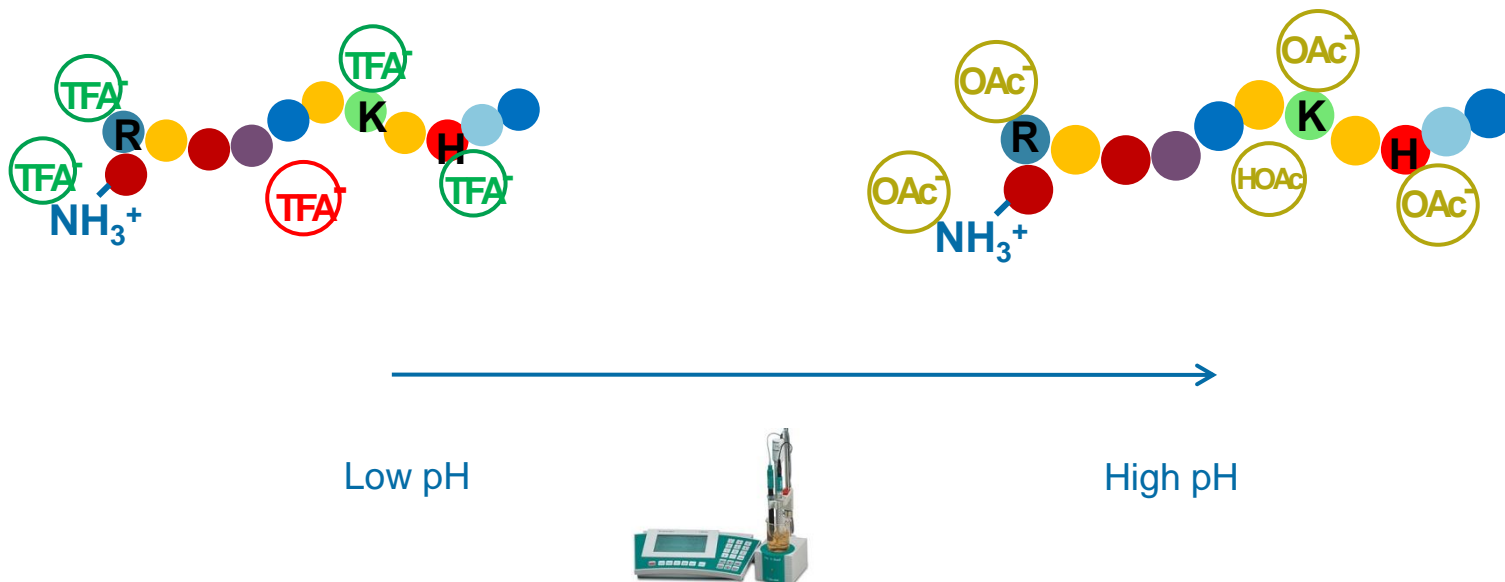
Solvent	Results (Dissolved or Undissolved)	Gross Peptide Concentration
Ultrapure water	Undissolved	N/A
PBS(pH 7.1 ± 0.1)	Dissolved	$\leq 1\text{mg/ml}$
DMSO	Dissolved	$\leq 10\text{mg/ml}$
Others*	N/A	N/A

pH Testing



◆ Why different pH value

- Free TFA acid may be present
- Different counter ions may result in various pH values of peptide solutions



Conclusions



Quantifying each component

Amino Acid Analysis
Peptide Content Analysis
Counter Ion Quantification Analysis
Moisture Content Analysis

Removing toxins

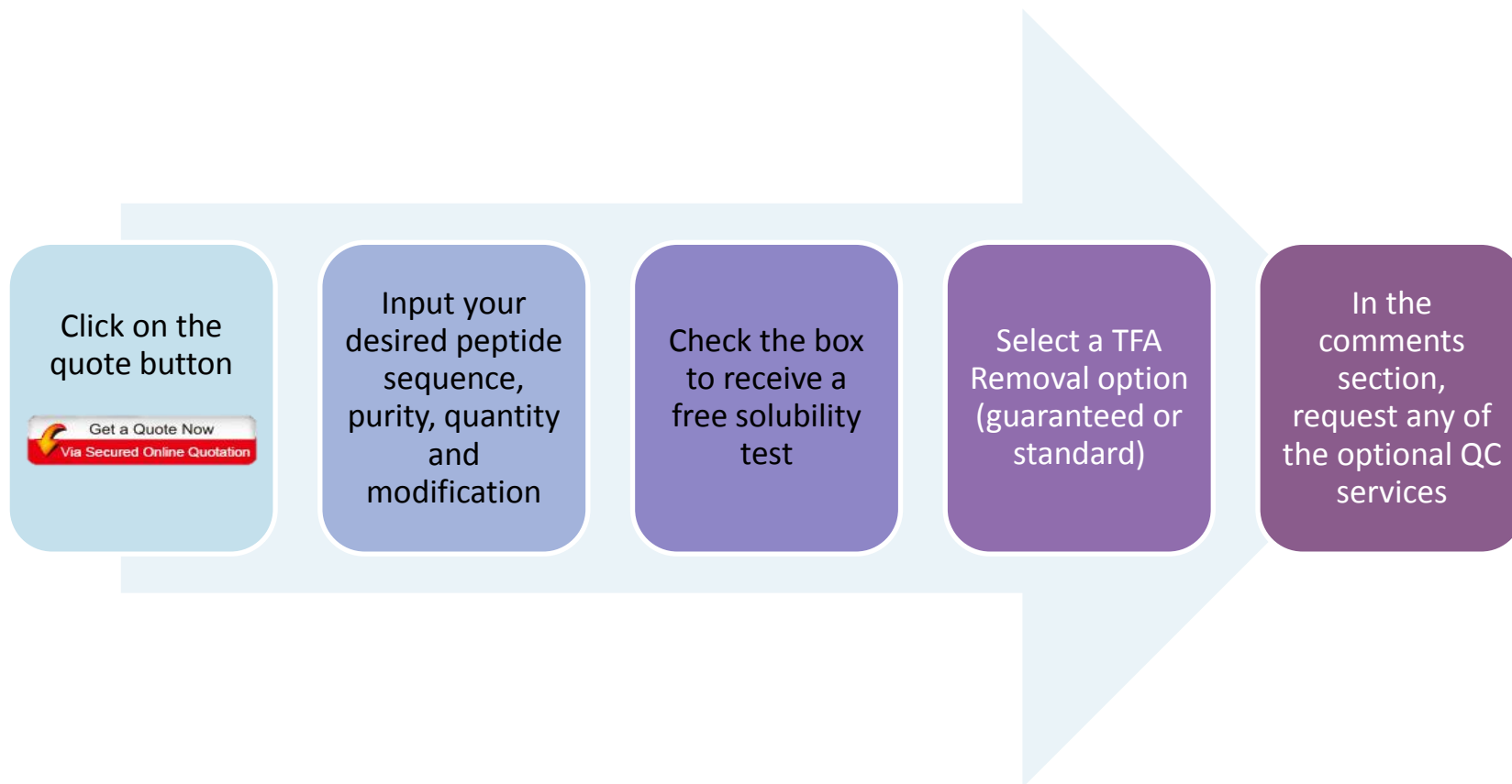
TFA Removal and Analysis
Endotoxin Analysis



Making sample prep easier

Solubility Test
pH Test

How can I request these services?



To learn more, visit

www.genscript.com/accuprep_quality.html

Peptide services at GenScript



Standard peptide synthesis

- Starting at \$3.2/AA
- Up to 200 AA
- Mg to kg quantity

Peptide Library

- Standard and micro-scale quantity
- Flexible purity options
- Peptide pooling available

Peptoid Synthesis

- Proteolytic resistant peptidomimetics
- Cost-effective, fast turnaround

Cosmetic Peptide Synthesis

- High batch-to-batch reproducibility
- High capacity

Click Peptide Synthesis

- O-acyl bond incorporation technology
- Increased peptide stability



www.genscript.com/peptide-services.html

*Thank you for your participation
We wish you success with your research*



Register for other webinars in the GenScript Webinar Series or download past webinars at <http://www.genscript.com/webinars.html>



Large scale genome editing for metabolic engineering of *E. coli*–
Yifan Li, Ph.D.

November 5, 2015, 9:00 AM or 2:00 PM EST



Optimizing soluble protein expression: codon optimization, RBS
design and expression vector design– *Rachel Speer, Ph.D.*

November 11, 2015, 9:00 AM or 2:00 PM EST



Antibody Drug Development: challenges & solutions – *Liusong Yin, Ph.D.*

November 18, 2015, 9:00 AM EST



If you have any other questions, visit www.genscript.com/faq_for_peptide
Or email: Lauren.Lu@genscript.com or laura.geuss@genscript.com