

# Identify the Optimal Purification Strategies for Your Recombinant Protein Production

Liyan Pang, Ph.D.

[liyan.pang@genscript.com](mailto:liyan.pang@genscript.com)



# Table of Contents

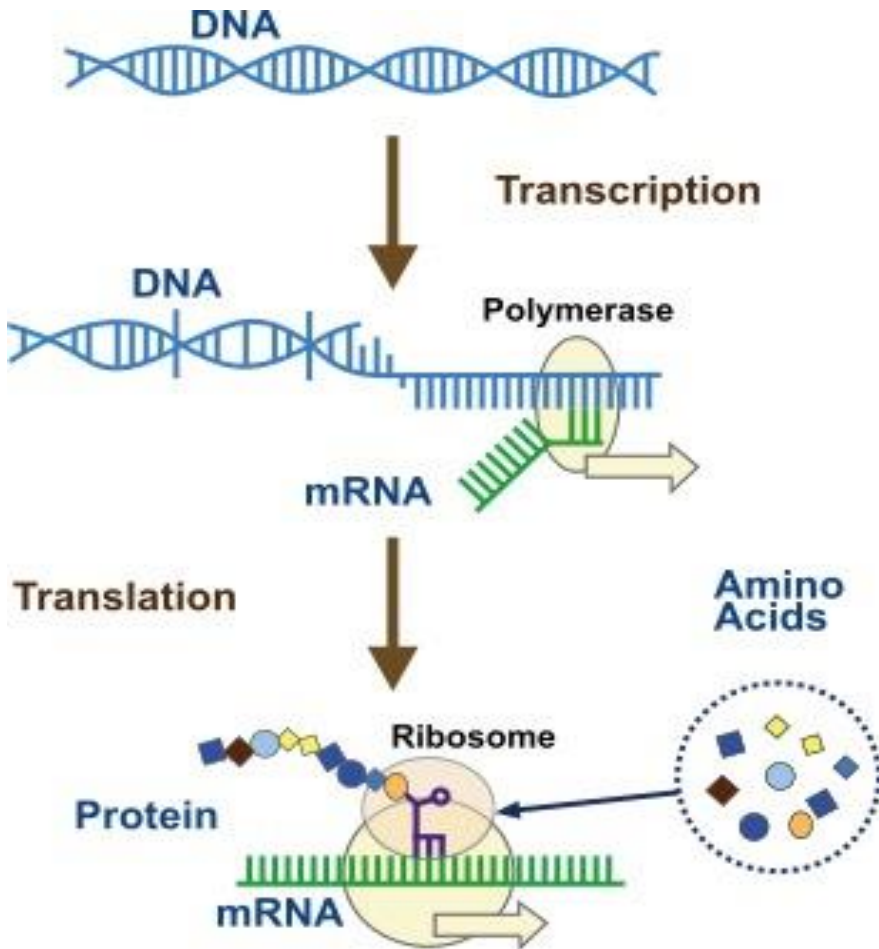


- 1 Protein Production Overview
- 2 Optimize Protein Expression
- 3 Protein Purification Strategies
- 4 Protein Tags
- 5 Endotoxin Removal
- 6 Advantages of GenScript Protein Services

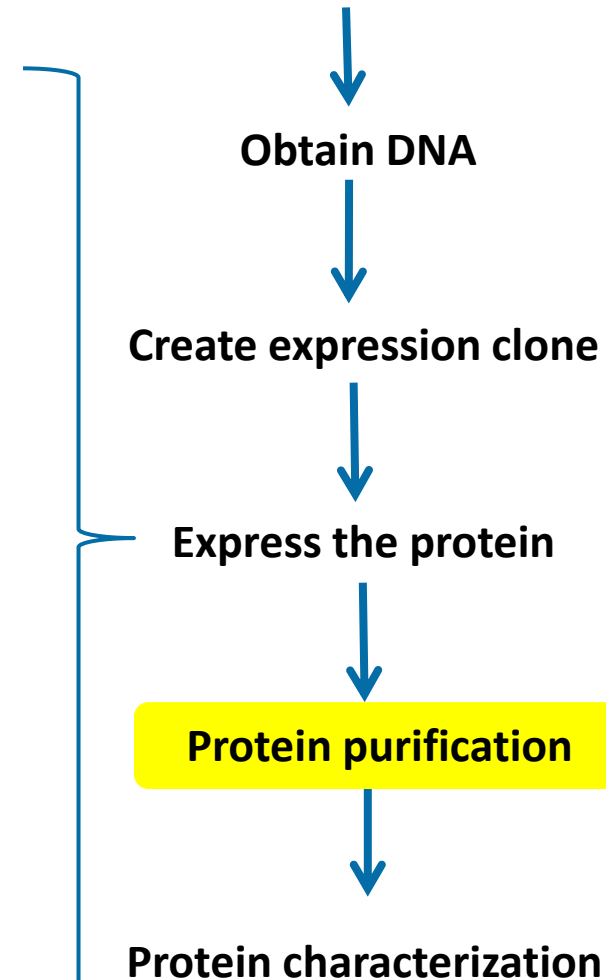
# Protein Production Workflow



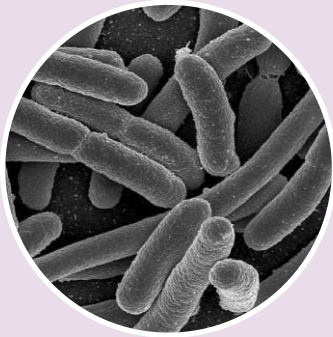
## Protein expression *in vivo*



## Protein name



# Expression Systems



## Bacteria

*E. coli*

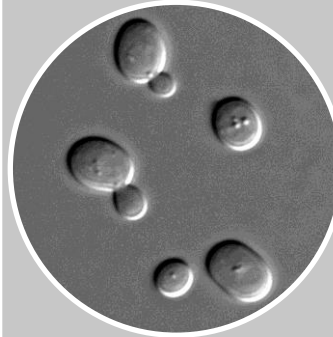
1. "Work horse"
2. Well established
3. High expression
4. Simple genetics
5. Easy scale up
6. Speed
7. Costs
8. Equipment



## Insect

*Sf9, Sf21, S2, High-5*

1. PTMs
2. Soluble proteins
3. High expressers



## Yeast

*S. cerevisiae*

*P. pastoris*

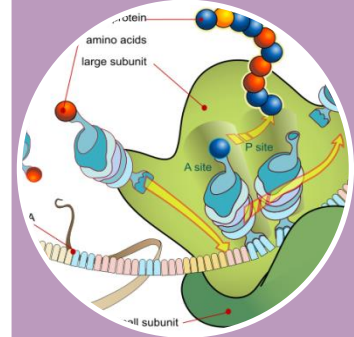
1. PTMs
2. Soluble proteins
3. High expresser



## Mammalian

*CHO, HEK, COS*

1. PTMs
2. Soluble proteins
3. Low expresser
4. Expensive



## Cell Free

*In vitro*

1. Expensive
2. Not reproducible
3. Scalability issues

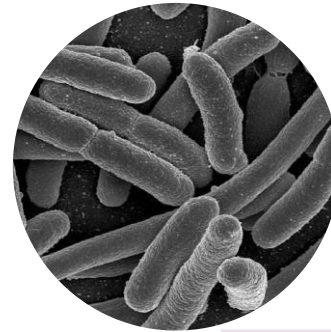


# Which Expression System to Choose?

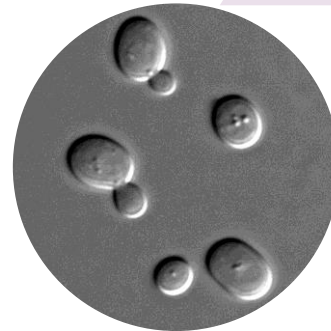


## Factors to Consider:

- ◆ Protein property
  - MW
  - Disulfide bonds
  - Post-translational modifications
  - Homogeneity
- ◆ Intended applications
  - Structural biology
  - Functional assays
  - Therapeutic protein/vaccines
  - Antigens for Ab production
  - Protein-protein interactions
- ◆ Yield
- ◆ Cost



Which to Choose



# GenScript Protein Production Services



From sequence to purified protein - gene synthesis included!

Expression System	Deliverables	Timeline	Price
<u><a href="#">BacPower™</a></u>	3mg purified protein guaranteed	6 -8 weeks	Starting from \$2,200
<u><a href="#">InsectPower™</a></u>	1mg purified protein guaranteed	8 -10 weeks	Starting from \$3,950
<u><a href="#">MamPower™</a></u>	3mg purified recombinant protein or 50mg purified antibody guaranteed	8 -12 weeks	Starting from \$8,499
<u><a href="#">YeastHigh™</a></u>	Customizable production up to 2000L	8 -10 weeks	Quote



## Eliminate the guesswork from your protein production work

- ◆ Evaluate whether your target protein expresses in your chosen system
- ◆ Identify the best expression system for your target protein
- ◆ Optimize your protein expression

Name	Service Type	Expression System(s)	Price	Timeline
<b>Standard</b>	Protein Expression Evaluation	<i>E. coli</i> (SC1653-B)	Starting from \$280	1-2 weeks
		Insect (SC1653-I)	Starting from \$400	3-4 weeks
		Mammalian (SC1653-M)	Starting from \$500	3-4 weeks
		All 3 Systems (SC1653-3S)	Starting from \$950	3-4 weeks
<b>Silver</b>	Protein Expression Optimization	<i>E. coli</i> (SC1667)	Login to inquire	2-3 weeks
<b>Gold</b>	Protein Expression Optimization	<i>E. coli</i> (SC1668)	Login to inquire	4-8 weeks

*GenScript's Solution for Expression Optimization*

# Case Study- Protein Expression Optimization



## ◆ Challenges:

- Pilot purification (final yield 1mg/L )- ~ 28kDa protein;
- Large amount of protein with large scale fermentation (1000L) and purification is needed.

## ◆ Strategies:

1. Expression improvement:
  - a. Promoter optimization
  - b. Strain optimization
  - c. pH optimization
  - d. Temperature and induction optimization
  - e. Inoculated quantity optimization
2. Recovery rate improvement during purification - Purification condition optimization
3. High-throughput screening with 48 conditions



# Case Study- Protein Expression Optimization



- ◆ 12X as much protein expression yield as original

1 mg/L starting protocol

2 mg/L T7 promoter/induction condition optimization

5 mg/L phoA promoter/induction condition optimization

6 mg/L growth condition optimization (pH)

12 mg/L seeding density optimization



- 1 Protein Production Overview
- 2 Optimize Protein Expression
- 3 Protein Purification Strategies
- 4 Protein Tags
- 5 Endotoxin Removal
- 6 Advantages of GenScript Protein Services

# What to Consider for Protein Purification?



**Goal:** To apply the optimal protein purification strategy specific to your recombinant protein, with which the highest level of purification is reached in the fewest steps.

- ◆ Protein characteristics
- ◆ Soluble vs. inclusion bodies
- ◆ Protein purification methods
- ◆ Protein with specific ligands
- ◆ Tagged protein
- ◆ Easy-to-degrade protein
- ◆ Concentration - Ultrafiltration
- ◆ Endotoxin removal

# Protein Expression as Inclusion Bodies



## ◆ What is an inclusion body?

- When *E. coli* is transformed to manufacture large amounts of recombinant protein, the protein sometimes forms dense aggregates of insoluble misfolded proteins, known as inclusion bodies.

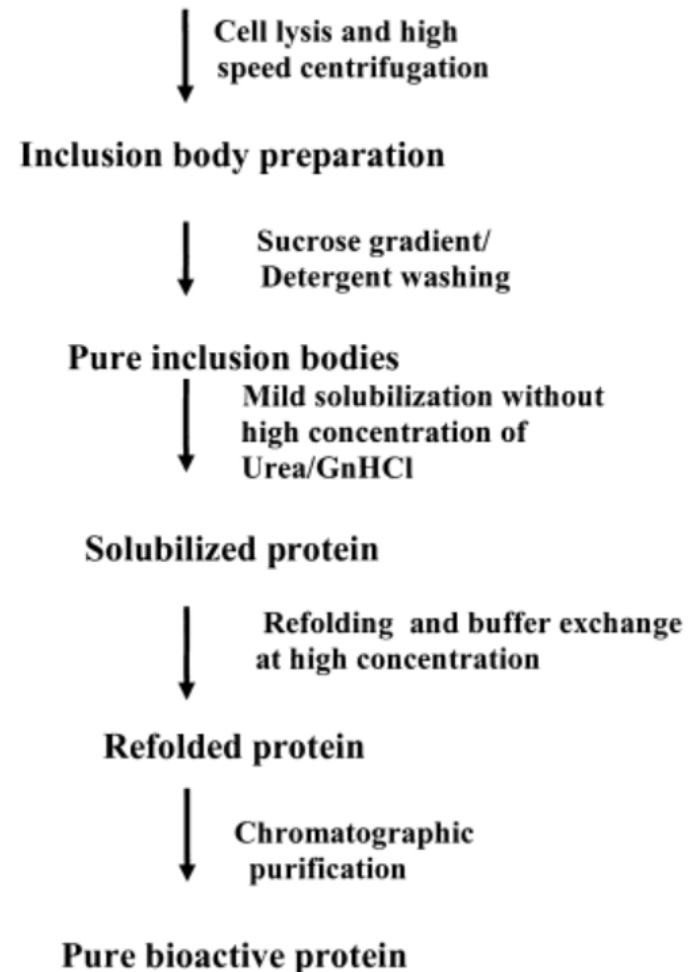
## ◆ Benefit

- allow high protein concentrations
- protect sensitive proteins from proteolytic (enzymatic) degradation
- protect the cell from any toxic proteins

## ◆ Challenge

- to solubilise and refold this protein into its correct 'active' form

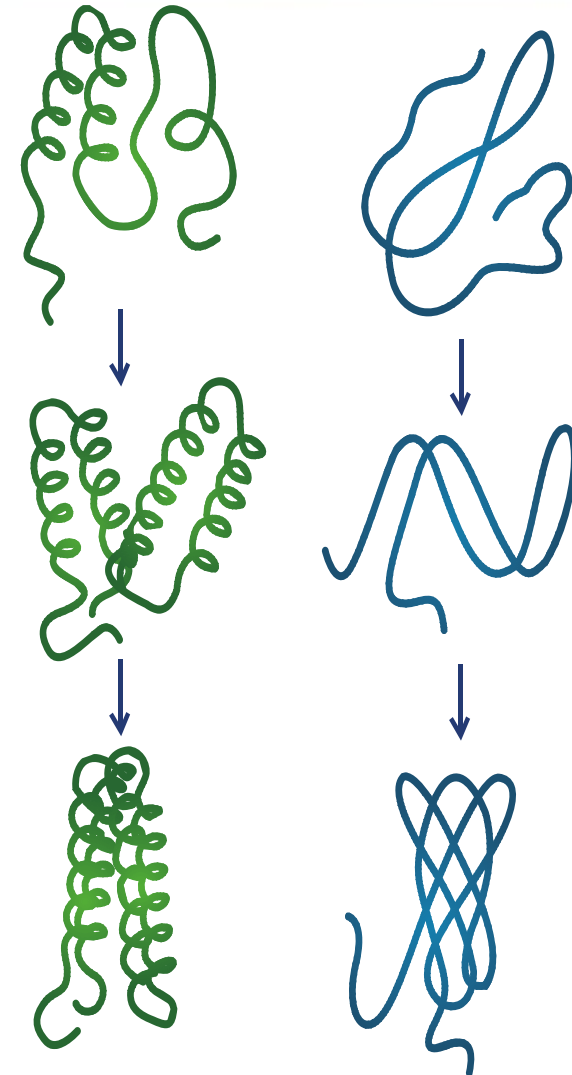
## *E. coli* expressing protein as inclusion bodies



# Protein Refolding Challenges



- ◆ All the information necessary for folding the peptide chain into its native structure is contained in the primary amino acid sequence of the peptide.
- ◆ There are vastly too many different possible conformations for a protein to fold by a random search.
- ◆ A new view of protein folding suggested that there is no single route, but a large ensemble of structures follow a many dimensional funnel to its native structure.
- ◆ Refolding conditions must be optimized for each individual protein.
- ◆ Important variables are:
  - buffer type
  - pH
  - ionic strength
  - Additives, often in combination (glycerol, redox reagents, saccharides, amino acids, metal ion, detergents, chaperones)



# GenScript's FoldArt™ Technology Overview



- Evaluation of target proteins' biochemical and biophysical properties
- Refolding optimizations
  - Selection of particular refolding strategy based on protein's sequence and the structural properties.
  - Buffer screening: Solutions for the inclusion body will be diluted to 20 different refolding buffers to determine which parameters affect the refolding results.
- Denaturant removal

Techniques: dilution, dialysis, diafiltration, gel filtration, and chromatography (ion exchange, size exclusion, and affinity)
- Validation

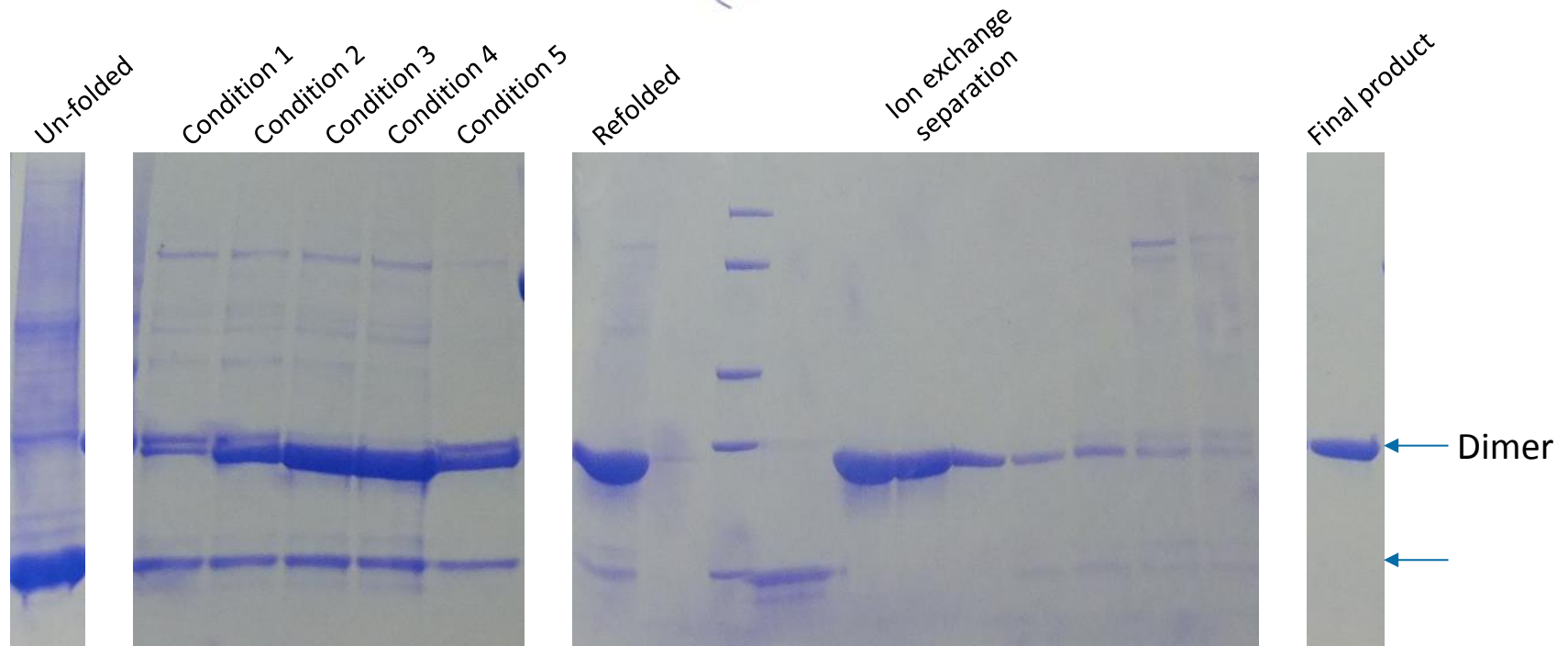
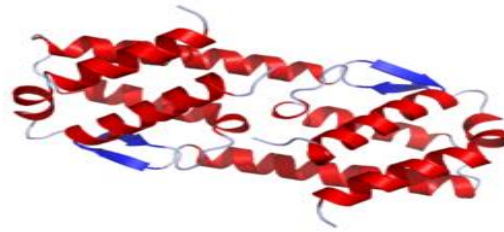
Refolding results will be validated by SDS-PAGE, HPLC and/or functional assay.

Over 95% of the inclusion bodies can be solubilized and refolded by our proprietary FoldArt™ protein refolding technology

# Case Study: Protein Refolding



- ◆ Human interleukin – 5: disulfide-bond linked homodimer as active form



# Protein Purification Methods



- Multiple choices on purification methods
  - Affinity column (GST, Ni-NTA, protein A/G/L resins, etc.)
  - Ion exchange (IEC)
  - Size exclusion (SEC)
  - Hydrophobic interaction chromatography (HIC)
- Double Tag strategy for big protein isolation





# Select Optimal Protein Purification Method



## ◆ Protein with special ligands

- IgG: Protein A or Protein G column;
- Fab & ScFv: Protein G or Protein L column;
- VhH/ScAb: Protein A column.

## ◆ Protein with tags

- Small tags:
  - His tag: Nickel-NTA column;
  - FLAG tag (DYKDDDDK): Anti-DYKDDDK conjugated resin;
- Big tags:
  - GST tag: Beads coated with Glutathione;
  - MBP tag: Amylose resin

## ◆ Tag free protein or additional purification

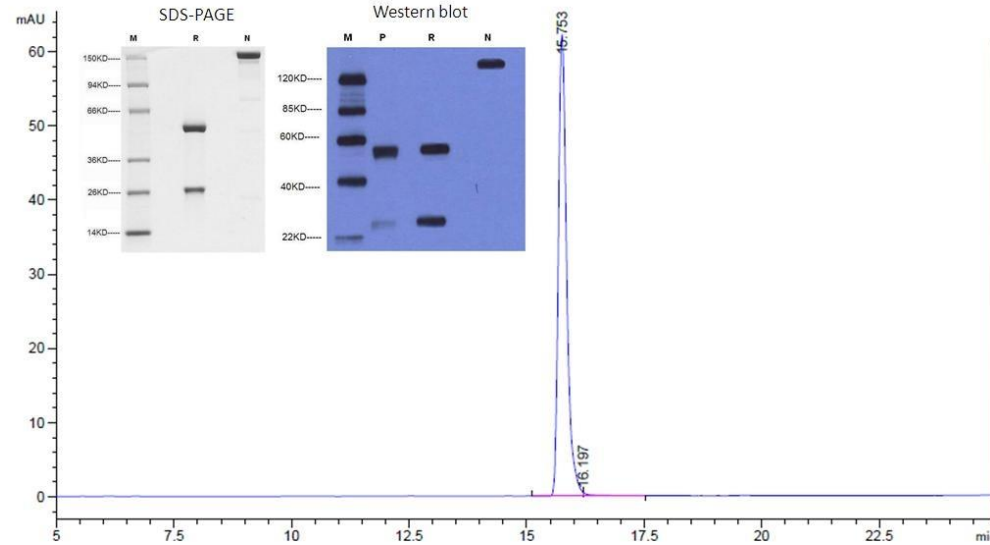
- Tag free protein: first to precipitate out the protein by ammonia sulfate.
- MW: proteins with distinct sizes can be separated by SEC.
- pI: IEC can be used to separate proteins based on their charges under particular pH.
- Protein with particular binding domain: Heparin column can be used to separate proteins with positive charged domains;
- Hydrophobicity: last resort to try, separation based on the proteins' hydrophobicity.

# Case Study: Recombinant mAb Production

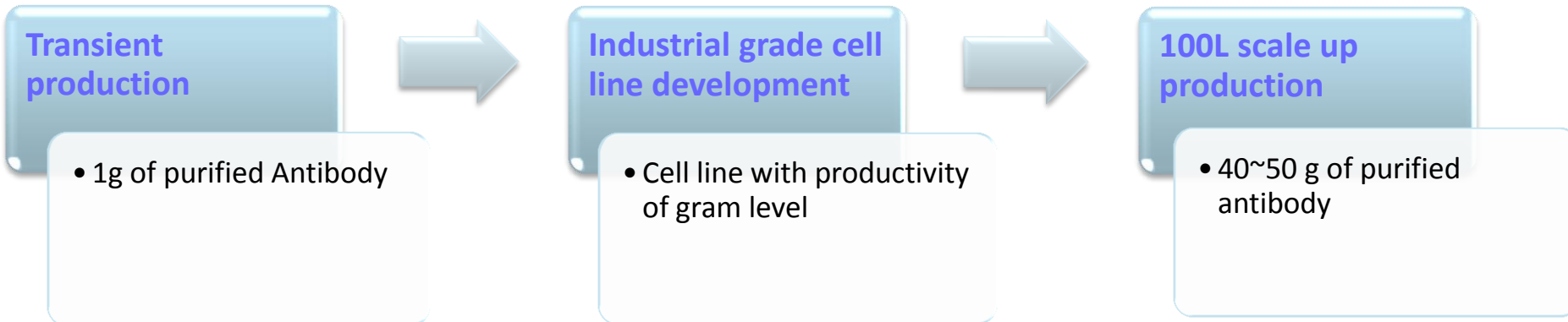


## Scale-up Production of a Recombinant mAb (hIgG1/kappa)

- Quality Standards:
  - Purity over 95% by SDS-PAGE and HPLC, even up to 99% purity
  - Aggregation less than 5%
  - Endotoxin Level: 0.05-1 EU/mg
  - Concentration: 20 mg/ml
  - Host cell protein 100 p.p.m.
  - Host cell DNA: < 100 pg/dosage



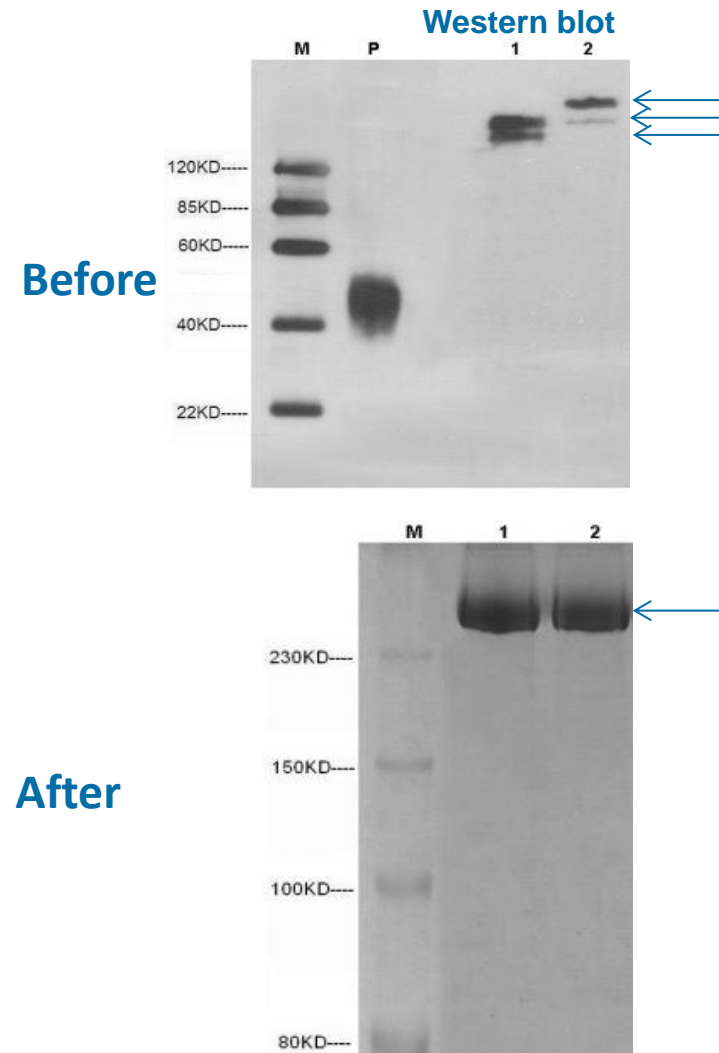
SDS-PAGE, Western blot and HPLC analysis



# Case Study: Prone-to-Degrade Protein



- ◆ Inconsistency of measured concentration
- ◆ Trouble shooting:
  - Transfection methods
  - Cell lysis
  - Purification
- ◆ Challenge: DNase is only partially responsible for the protein degradation. This protein itself is prone-to-degrade.
- ◆ Solutions:
  - Remove DNase
  - Add protease inhibitor to every step
  - Optimize buffer components
  - Add protein stabilizers
  - Lyophilization immediately after protein purification
  - Storage temperature

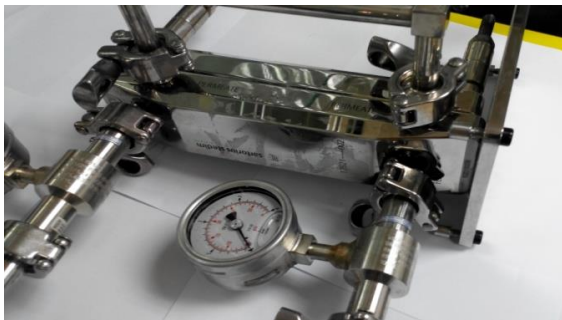


# Protein Concentration - Ultrafiltration



## Cross-flow filtration

### Plate-and-Frame Module



### Advantages

- High allowable work pressure
- Easy to clean
- Easy to replace membranes

### Hollow-Fiber Module



## Dead-end filtration



### Advantages

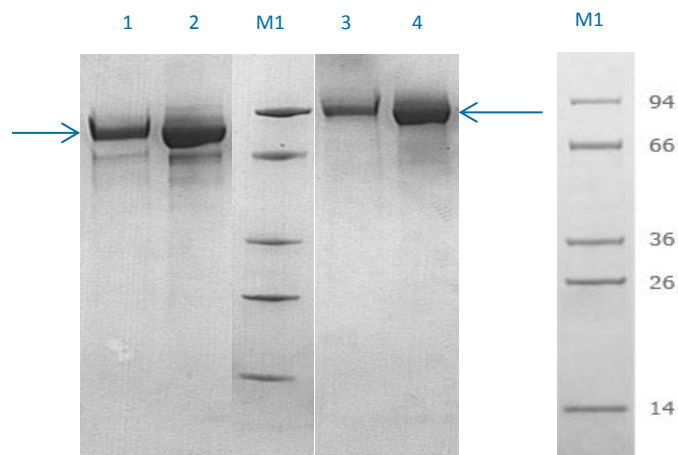
- Ultra-fast spin times (process 15mL samples as fast as 10-15 minutes)
- Ultra-high recoveries (typically >90%)
- Relative low costs

# Case Study on Protein Concentration



Slice Cassettes; MWCO:30 KD; Membrane Area:0.1m<sup>2</sup>

Sample	MW (KD)	Initial concentration (mg/ml)	Initial volume (L)	Final concentration (mg/ml)	Final volume (ml)	Tem (°C)	Recovery (%)
Protein S	79.5	0.05	5	1.000	243	RT	97.2
Protein Q	92.7	0.04	5	0.800	196	RT	98.0



M1: Protein Marker

Lane1: The **protein S** before Concentration (30ul)

Lane2: Concentrated **protein S** (3ul)

Lane3: The **protein Q** before Concentration (30ul)

Lane4: Concentrated **protein Q** (3ul)



- 1 Protein Production Overview
- 2 Optimize Protein Expression
- 3 Protein Purification Strategies
- 4 Protein Tags
- 5 Endotoxin Removal
- 6 Advantages of GenScript Protein Services



## ◆ Protein information:

- Fusion protein (PI, Molecular weight, concentration)
- Tag type
- Protease cleavage site
- Tag-free protein (PI, Molecular weight)

## ◆ Protease information:

- Protease activity
- Protease reaction condition

# Affinity Chromatography with Tagged Proteins



## ◆ Affinity tags

- His
- GST
- MBP
- Flag
- SUMO

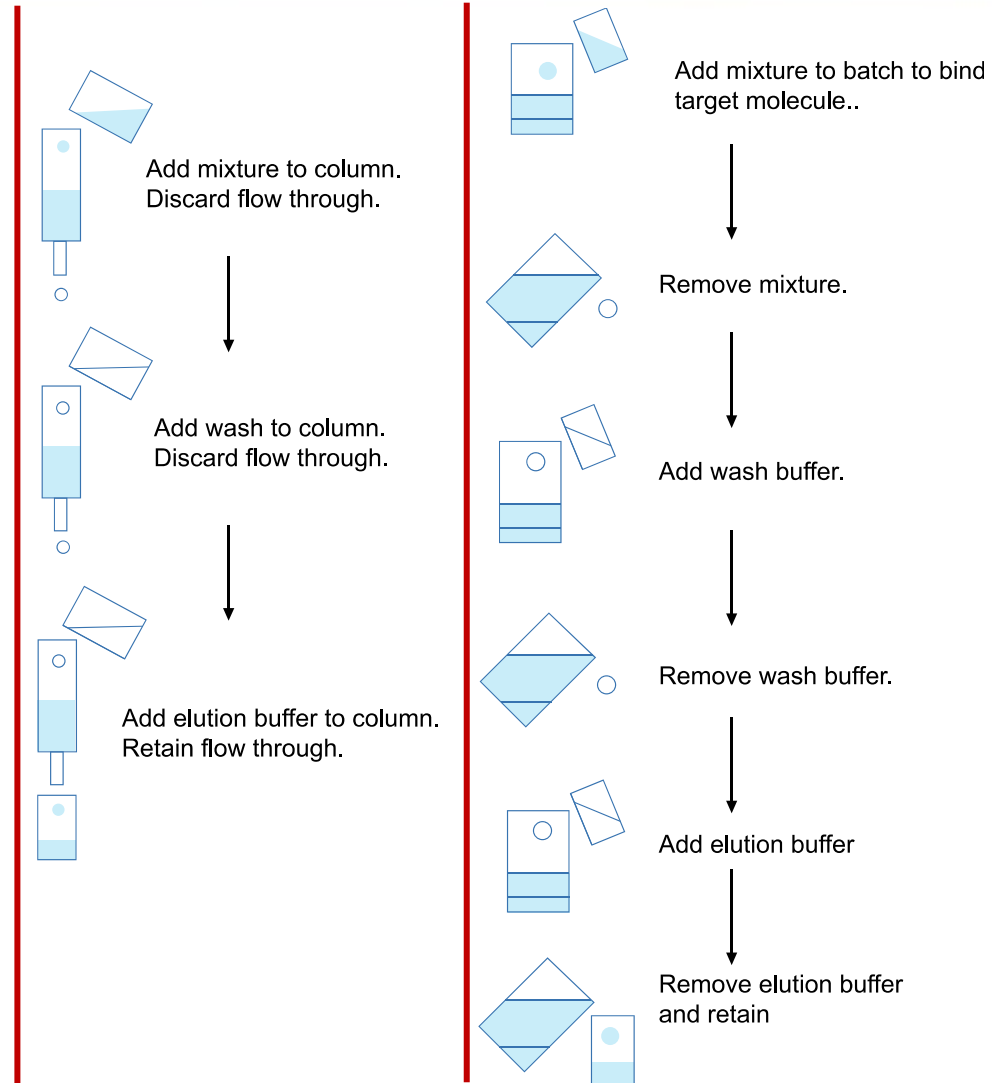
## ◆ Matrices & elution conditions

## ◆ Methods

- Column
- Batch

## ◆ Tag removal by site-specific protease

- TEV
- Enterokinase
- SUMO protease



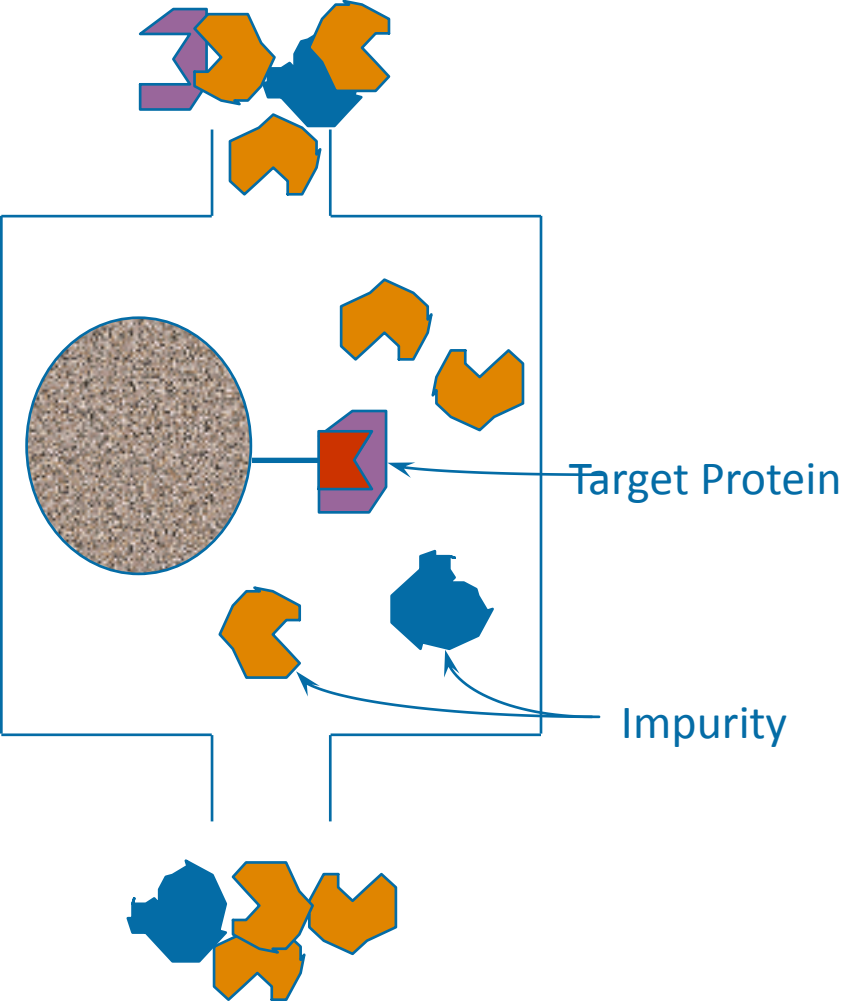
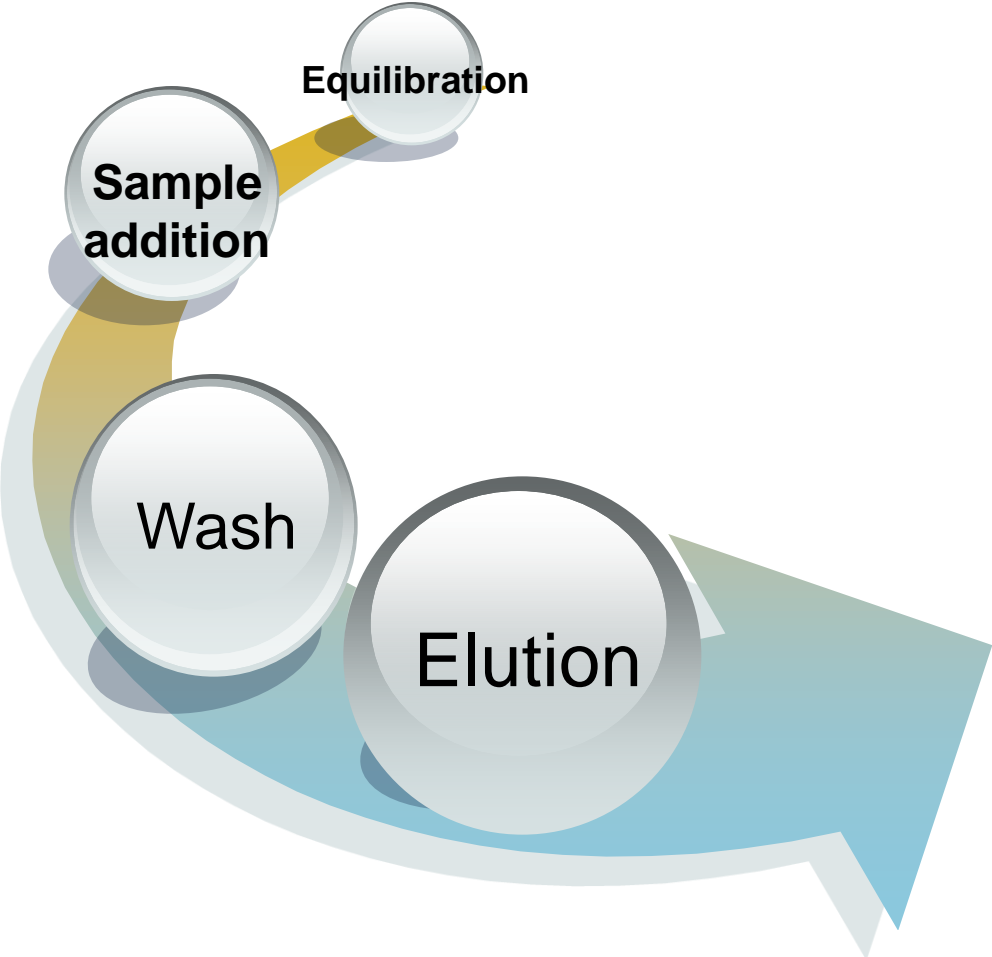


# Commonly Used Protein Tags

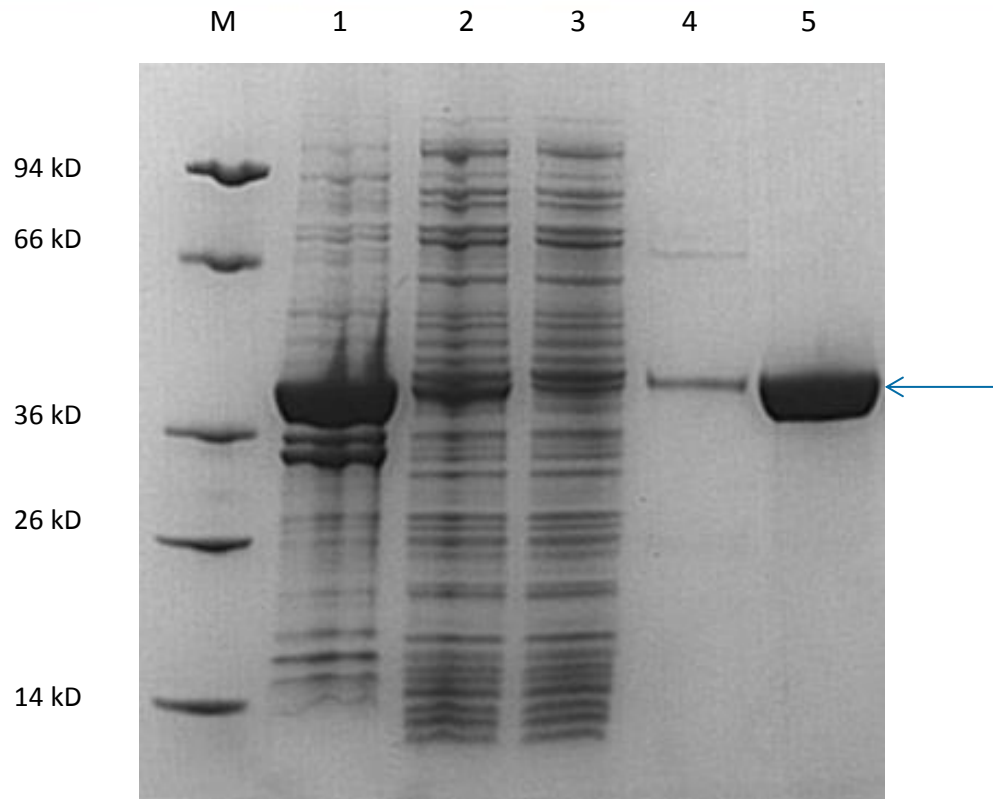


Tags	Amino Acids	Size (KDa)	Matrix	Comments
His	6-8	~1	Nickel	Most commonly used
FLAG	8	~1	Anti-FLAG	Specific
MBP	396	40	Amylose	↑ solubility; periplasm
GST	211	26	Glutathione	Dimerization
NusA	495	54.9	N/A	↑ solubility
Trx	109	11.7	N/A	↑ Disulfide bond
DsbA	208	23.1	N/A	↑ Disulfide bond; periplasma
SUMO	101	11.6	N/A	↑ solubility
Z-tag	91	10.6	Protein A	↑ solubility

# Workflow of Tagged Protein Purification



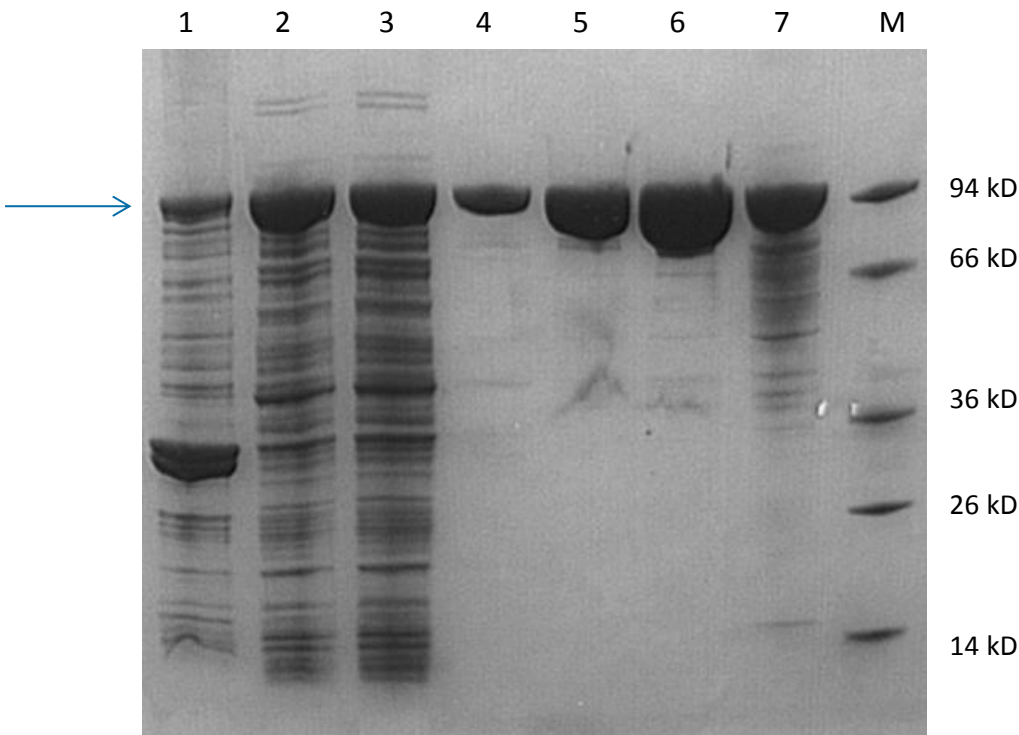
# Case Study: One Step His-tag Nickel Purification



Lane:

- 1: Precipitation after cell lysate centrifugation
- 2: Supernatant after cell lysate centrifugation
- 3: Flow through
- 4: Wash with 50 mM imidazole
- 5: Elute with 500 mM imidazole

# Case Study: One Step MBP-Amylose Purification



- Lane:
- 1: Precipitation after cell lysate centrifugation
  - 2: Supernatant after cell lysate centrifugation
  - 3: Flow through
  - 4-6: Elute with amylose gradient
  - 7: Resin

# Commonly Used Proteases

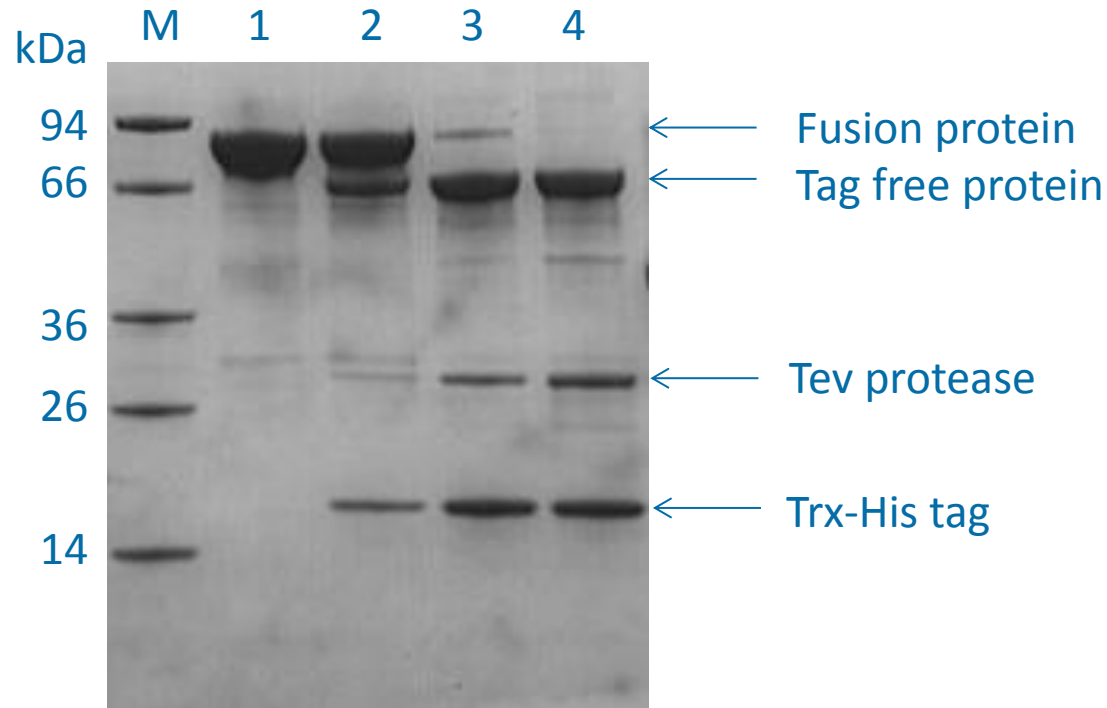


Protease	Recognition sequence	Amino acid residues	Temperature range (°C)	pH Range	Molecular weight
Tev	ENLYFQ <sup>^</sup> G	G	4~30	5.5~8.5	28KDa
EK	DDDDK <sup>^</sup>	None	4~37	6~9	22.7KDa
Thrombin	LVPR <sup>^</sup> GS	GS	4~37	~7	36KDa
3C	LEVLFQ <sup>^</sup> GP	GP	4~15	7~8.5	42KDa
SUMO	Small ubiquitin-like modifier	None	4~30	5.5~9.5	25KDa,28KDa

# Case Study: Trx Tag Removal by TEV Protease



4°C, overnight



SDS-PAGE analysis of Tev protease digestion of Trx-His protein followed by Coomassie Blue Staining



- 1 Protein Production Overview
- 2 Optimize Protein Expression
- 3 Protein Purification Strategies
- 4 Protein Tags
- 5 Endotoxin Removal
- 6 Advantages of GenScript Protein Services

# Endotoxin Removal



## ◆ What are endotoxins?

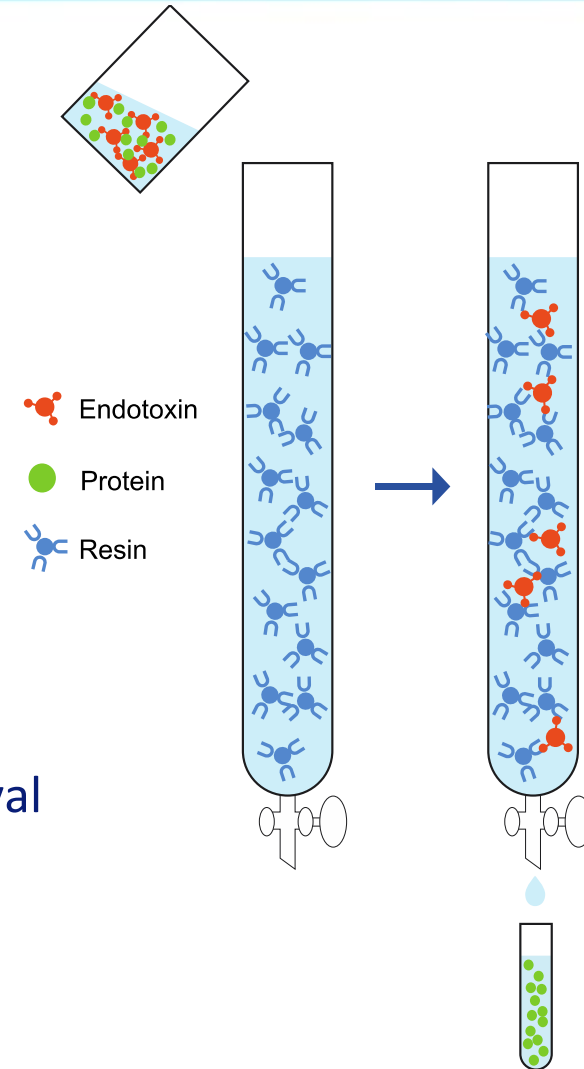
- Endotoxins, also known as lipopolysaccharides (LPS), are large molecules found in the outer membrane of Gram-negative bacteria, which elicit strong immune responses in animals.

## ◆ Detection

- Gel clot method
- Chromogenic method

## ◆ Removal methods:

- Polymyxin B (PMB) – affinity based ToxinEraser™ (L00338) by GenScript allows highly efficient removal of endotoxin down to 0.1 EU/ml
- Size exclusion chromatography (SEC)
- Ion exchange chromatography (IEC)







# Case Study: Endotoxin Removal by PMB

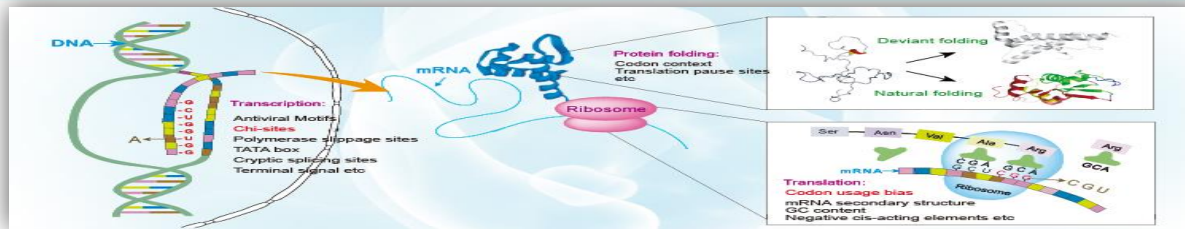
GenScript Endotoxin Removal Services: Endotoxin  $\leq 1$  EU/ $\mu$ g;  $\leq 0.1$  EU/ $\mu$ g;  $\leq 0.01$  EU/ $\mu$ g

	Protein 1	Protein 2	Protein 3
Volume of resins	3 ml	3 ml	3 ml
Volume of sample	15 ml	15 ml	15ml
Initial endotoxin	500,000 - 2,000,000 EU/ml	> 40,000,000 EU/ml	> 40,000,000 EU/ml
Final concentration	2.2 mg/ml	1 mg/ml	0.8 mg/ml
Final endotoxin	64 – 128 EU/ml	20 – 40 EU/ml	12.5 – 25 EU/ml
Final endotoxin	0.029 – 0.058 EU/ $\mu$ g	0.02 – 0.04 EU/ $\mu$ g	0.016 – 0.032 EU/ $\mu$ g



- 1 Protein Production Overview
- 2 Optimize Protein Expression
- 3 Protein Purification Strategies
- 4 Protein Tags
- 5 Endotoxin Removal
- 6 Advantages of GenScript Protein Services

# Advantages of GenScript Protein Services



## ◆ Core in-house technologies for expression optimization & production efficiency.

- OptimumGene™ – expression system specific codon optimization
- BacPower™ – increase bacterial soluble expression
- FoldArt™ – ensure appropriate protein refolding
- YeastHigh™ – high copy-number gene selection technique
- BacuVance™ – for protein secretion from baculovirus-infected insect cells
- MamPower™ – technology licensed from NRC for rapid recombinant protein production with high yield
- DoubleTag strategy – for big protein isolation

# Advantages of GenScript Protein Services



## ◆ One-stop service from sequence to purified proteins with large capacity.

### ▶ Guaranteed Protein Expression Package

- 3 mg purified soluble protein from \$2,200
- Subcloning
- Transformation
- Expression
- Refolding



### ▶ Protein Expression and Purification Services

- Bacteria
- Yeast
- Baculovirus/insect cells
- Mammalian

### ▶ Large-scale Protein Production Services

- Bacterial fermentation up to 1,000 L
- Yeast fermentation up to 500 L
- Baculovirus/insect cell production up to grams
- Mammalian cell production up to grams



### ▶ Bioprocessing Services

- Mammalian protein expression services
- Stable cell line development & protein production

**OptimumGene™ Gene Design Service**

— Minimum 10-fold increase in protein express level

# Advantages of GenScript Protein Services



- ◆ Flexible production scales
- ◆ Fast turn-around time (from sequence to purified protein in as little as 4 weeks)



## Capacity:

Bacteria	Yeast	Baculovirus	Mammalian
1,000 L	500 L	100 L	500 L
Fermentor	Fermentor	Wave™ Mixer	Wave™ Mixer Sartorius bioreactor Hyclone SUB bioreactor

# Variety of GenScript Protein Services



## 💧 Protein Expression & Purification

Bacteria, Insect, Mammalian, Yeast,  
Customized protein services, Fermentation,  
Transient expression, Stable cell lines

## 💧 Large Scale Production

Scale up production in Bacteria, Yeast, Insect  
and Mammalian cells, Large-scale  
Recombinant Mab Production

## 💧 Other Protein Services

Structural Biology, HTP Variants, Chemical  
Protein Synthesis, Endotoxin removal, Codon  
optimization, Custom protein purification,  
Protein characterization, Protein Refolding

## 💧 Expression Evaluation & Optimization

Small scale expression testing and optimization in  
bacterial, insect and mammalian expression  
systems

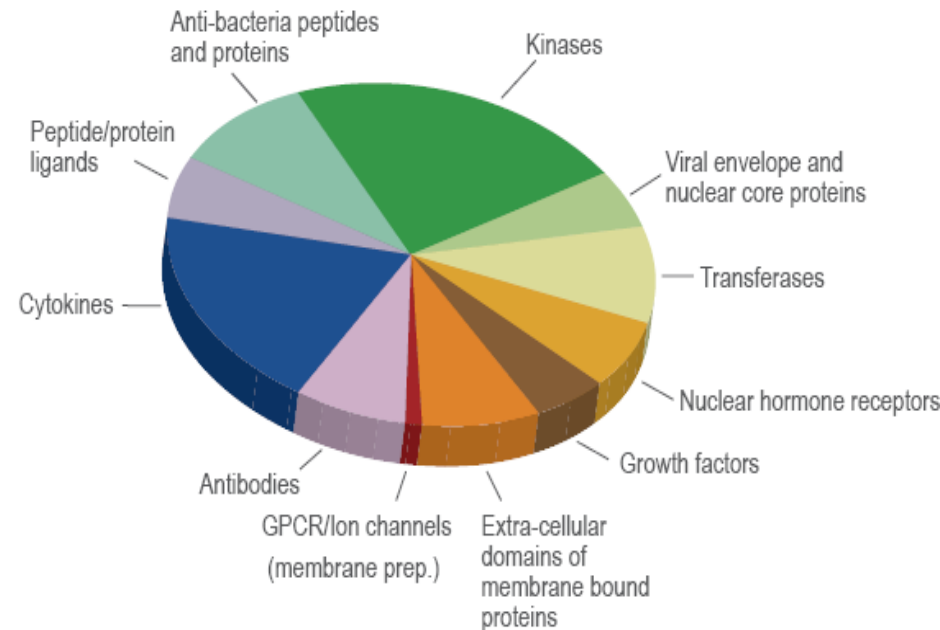
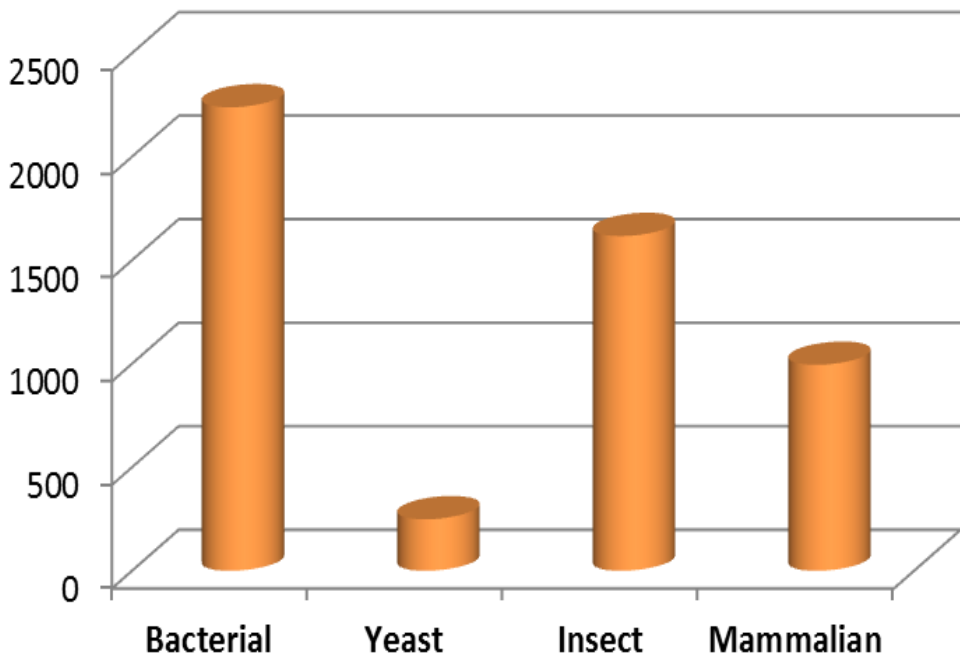
## 💧 BioProcess

Upstream & Downstream Process development,  
GLP-compliant Bioprocess Services

## 💧 Resources and Technical References

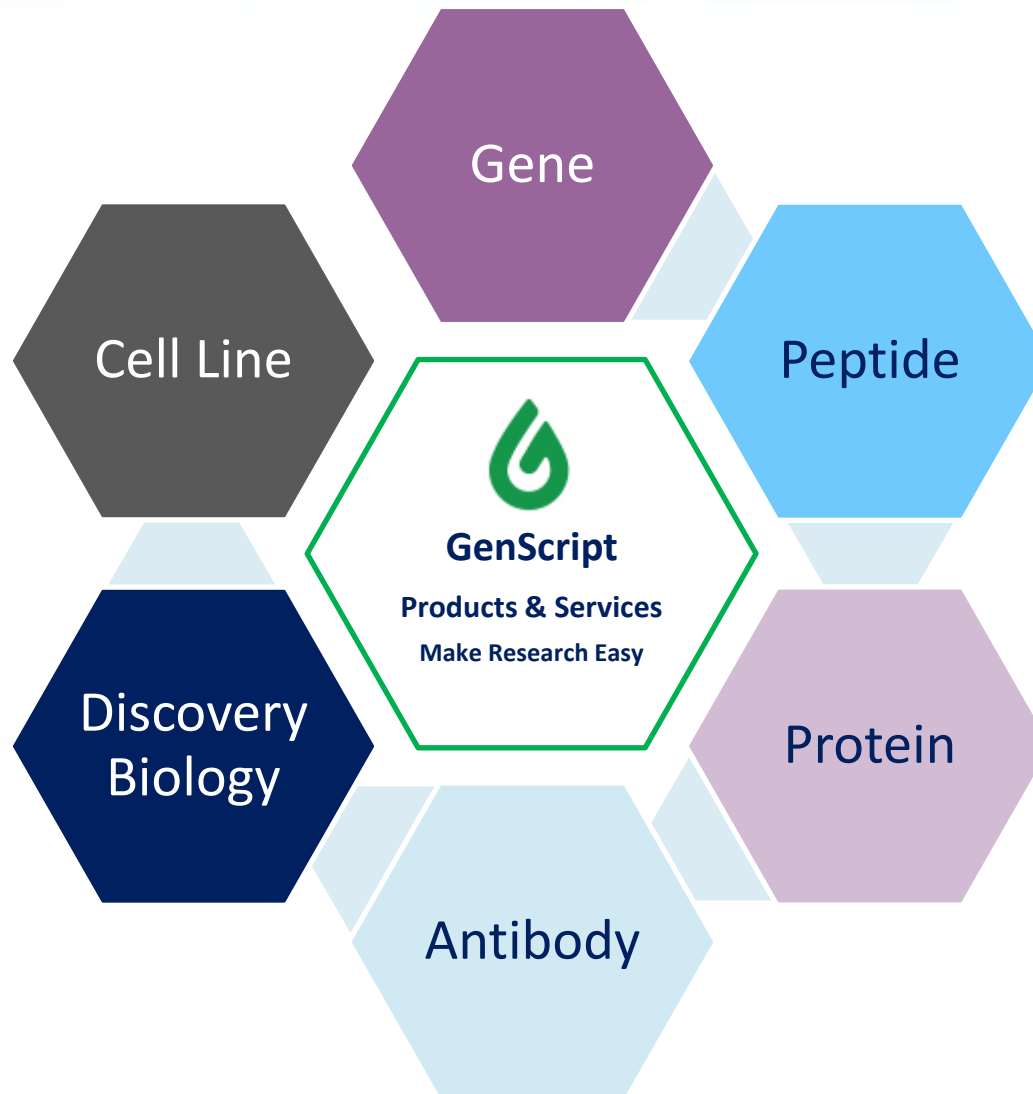
PSORT II, WoLF PSORT II, Case Studies,  
Downloads, FAQs, ProtBank™ Protein Database  
Protein News Page, Publications, Service  
Selection Guide

# GenScript's Experience in Protein Expression & Purification



GenScript has delivered over **5,000** proteins in four expression systems. Statistics showed **95%** success rate for all protein projects.

# About GenScript





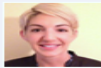
Thank you for your participation  
We wish you all success in your Research



Register for other webinars in the GenScript Webinar Series @  
<http://www.genscript.com/webinars.html>

#### Anti-idiotypic antibodies – A powerful tool for antibody drug studies

Need a reliable tool for your antibody drug PK/PD or immunogenicity studies? Anti-idiotypic antibodies (anti-IDs) serve as powerful tools for such studies, as they are able to differentiate antibody drug from endogenous antibodies and allow tracking of antibody drug in biological fluid. However, it has been challenging to develop highly specific and sensitive anti-IDs due to the low percentage of anti-IDs relative to other naturally-occurring antibodies in immunized animals. This webinar will take an in-depth look into the field of anti-IDs, including the obstacles and solutions surrounding the generation of highly sensitive, highly specific anti-IDs, as well as their use in vaccine development, PK/PD and immunogenicity studies.



Presented by: Michelle Parker, Ph.D.

November 20, 2014/  
8:00 am

Register now

November 20, 2014/  
2:00 pm

Register now

#### Peptide design strategy: basics, optimization, and application

Is your peptide sequence optimally designed for your unique research application? Peptide design effects the overall stability and usability of your peptide in experiments, as well as the turnaround time for your custom peptide. In this session, we will review the basics of peptide design strategy including sequence design, modification selection, and purity choice. Case studies, best practices, and design tools will be introduced. In addition, we will explore specific trends in peptide design as it applies to research applications including drug discovery, drug delivery, structural biology, and biomaterials.



Presented by: Tiffany Gupton Campolongo, Ph.D.

November 25, 2014/  
8:00 am

Register now

November 25, 2014/  
2:00 pm

Register now

#### Fusion partner for recombinant soluble protein production in *E. coli*

Are you working on an *E. coli* expression project and looking for ways to generate soluble protein? If so, this interactive webinar may be for you. Target protein solubility is one of the most important end goals of a recombinant protein production campaign yet it remains a significant bottleneck. While there are various ways to optimize expression in order to produce soluble protein, a fusion partner based approach can be very useful. This webinar reviews the practical aspects of a fusion partner strategy. Topics include but are not limited to - the need for protein solubility, the relationship between solubility and protein functionality, use of soluble fusion partners, examples of commonly used partners, identifying novel fusion partners and pros and cons of a fusion partner approach.



Presented by: Keshav Vasanthavada

December 3, 2014/  
8:00 am

Register now

December 3, 2014/  
2:00 pm

Register now