

Recombinant Protein Expression & Purification -- Challenges & Solutions

Liyan Pang, Ph.D.

liyan.pang@genscript.com



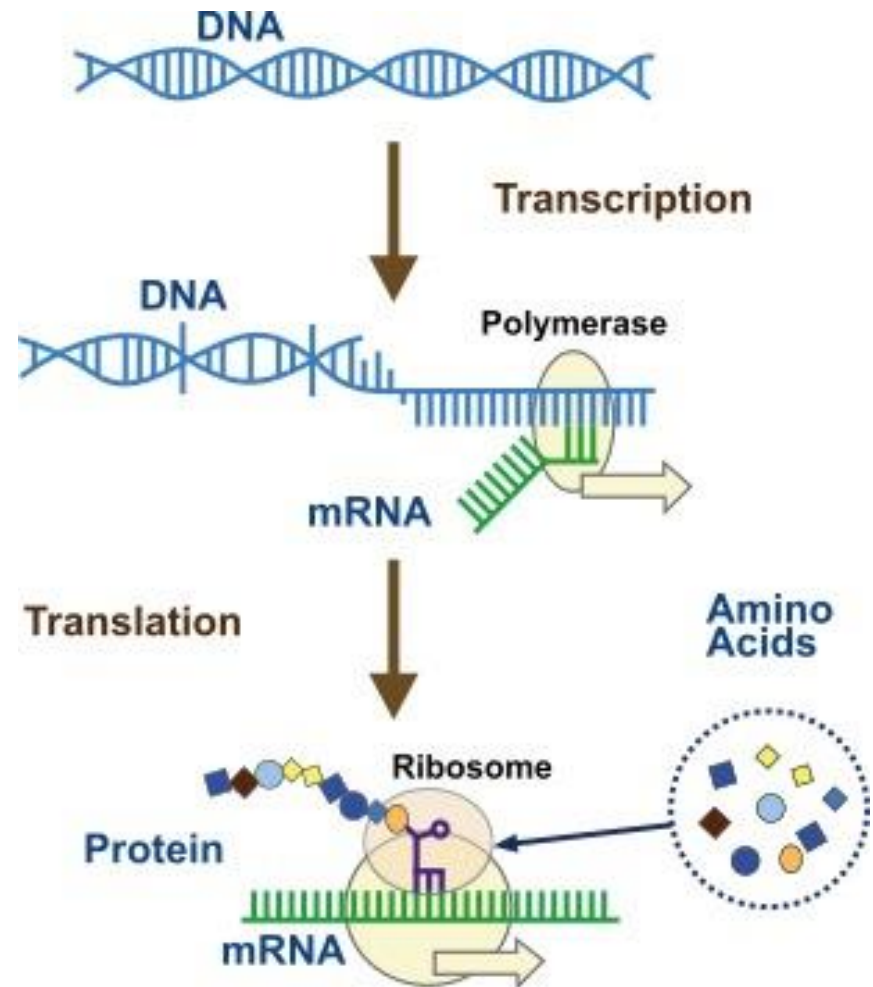


- 1 Choose Expression System
- 2 Optimize Protein Expression
- 3 Protein Refolding
- 4 Protein Purification
- 5 Difficult-to-Express Proteins
- 6 Advantages of GenScript Protein Services

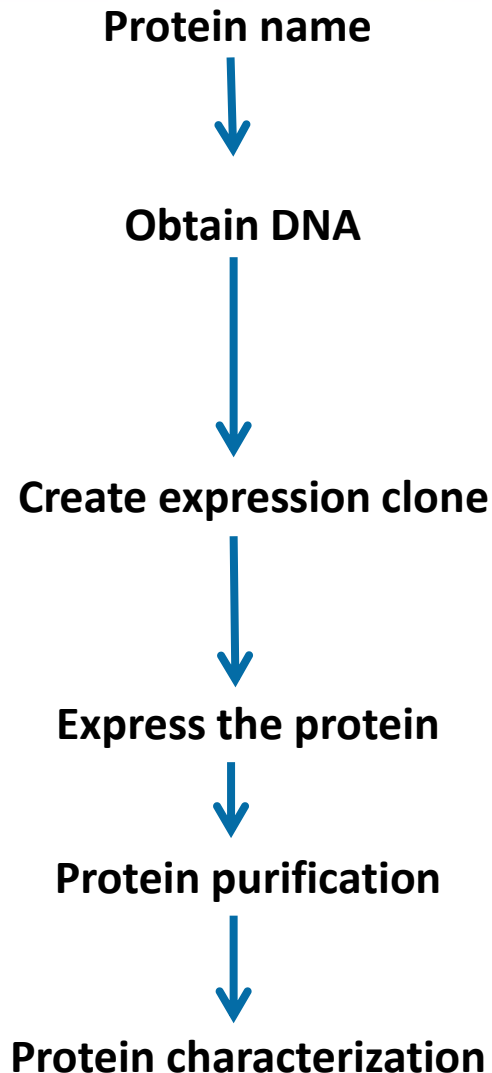
Challenges in Protein Expression



- ◆ Soluble expression
- ◆ Correct conformation
 - Disulfide bond
 - Chaperonins
 - Refolding
- ◆ Production efficiency
- ◆ Purity
- ◆ Difficult-to-express proteins
 - Membrane protein
 - Toxic protein
 - Easy-to-degrade protein

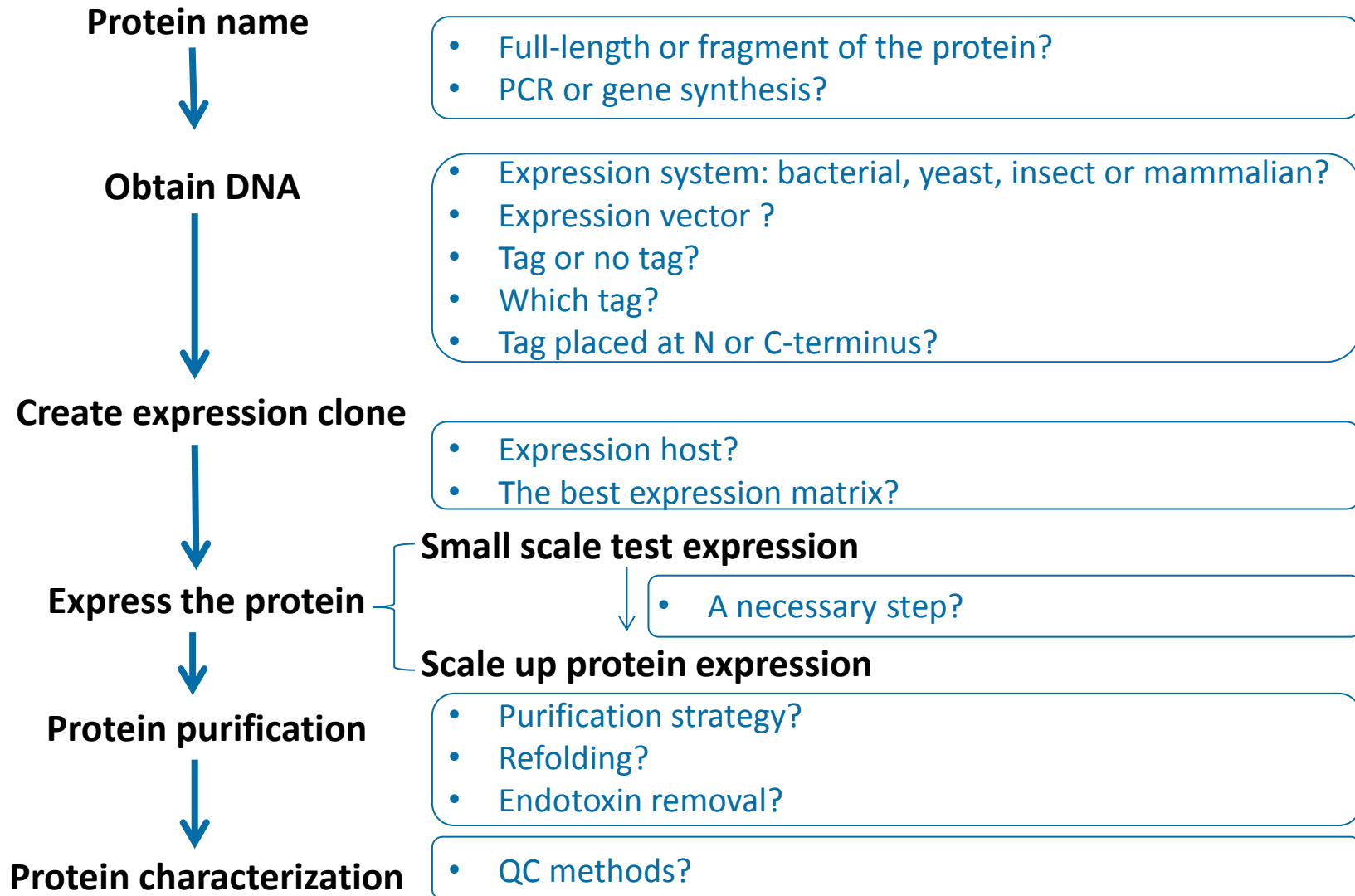


Protein Expression Impacting Factors



- ◆ Sequence
- ◆ Expression system
 - Bacterial
 - Yeast
 - Insect
 - Mammalian
- ◆ Vector
 - Promoter
 - Tag
- ◆ Host strain
- ◆ Expression conditions
 - Medium component
 - Temperature
 - Inducer concentration & time
 - Inoculation volume

Protein Expression Common Questions



Expression Systems



Bacteria

E. coli

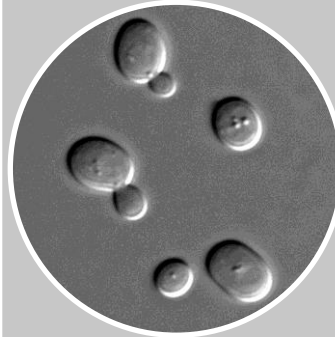
1. "Work horse"
2. Well established
3. High expression
4. Simple genetics
5. Easy scale up
6. Speed
7. Costs
8. Equipment



Insect

Sf9, Sf21, S2, High-5

1. PTMs
2. Soluble proteins
3. High expressers



Yeast

S. cerevisiae
P. pastoris

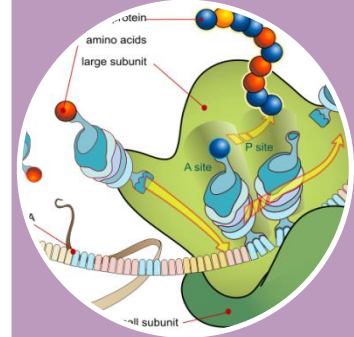
1. PTMs
2. Soluble proteins
3. High expresser



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

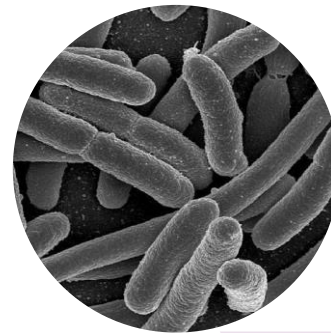


Which Expression System to Choose?

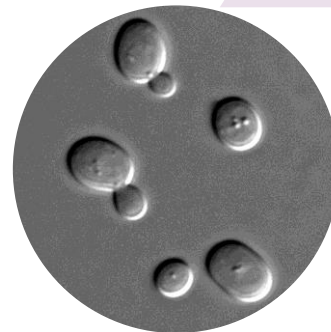


Factors to Consider:

- ◆ 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



Expression System Selection



Expression System	Pros	Cons	Intended Applications
<u>Bacterial</u>	<ul style="list-style-type: none"> Relatively inexpensive Simple genetics Easy to manipulate Easy scale up Fast expression 	<ul style="list-style-type: none"> Lack of efficient post-translational modifications Codon usage issues Inclusion bodies Low yield and activity for some eukaryotic proteins Difficult to express higher MW proteins 	<ul style="list-style-type: none"> <u>Structural biology</u> <u>Functional assays</u>
<u>Yeast</u>	<ul style="list-style-type: none"> Diverse post-translational modifications Low cost of culture media Industry-scale fermentation 	<ul style="list-style-type: none"> Improper glycosylation Excessive glycosylation 	<ul style="list-style-type: none"> <u>Antigen production</u>
<u>Insect</u>	<ul style="list-style-type: none"> Good secretion Post translational modifications resemble mammalian system Suitable for toxic gene products 	<ul style="list-style-type: none"> Long production time Relative high media costs 	<ul style="list-style-type: none"> Protein-protein interaction (Bacterial & Yeast expression systems recommended)
<u>Mammalian</u>	<ul style="list-style-type: none"> Comprehensive post-translational modifications Excellent method for the production of bioactive proteins 	<ul style="list-style-type: none"> Long production time High media costs Protein yields relatively lower 	<ul style="list-style-type: none"> Therapeutic protein (Yeast & Mammalian expression systems recommended)

GenScript Protein Standard Services



From sequence to purified protein - gene synthesis included!

Expression System	Deliverables	Timeline	Price
<u>BacPower™</u>	3mg purified protein guaranteed	6 -8 weeks	Starting from \$2,200
<u>InsectPower™</u>	1mg purified protein guaranteed	8 -10 weeks	Starting from \$3,950
<u>MamPower™</u>	3mg purified recombinant protein or 50mg purified antibody guaranteed	8 -12 weeks	Starting from \$8,499
<u>YeastHigh™</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
 - PROtential™ Standard packages
 - Before scale-up protein production, to avoid waste on your time & valuable resources
- ◆ Optimize your protein expression
 - PROtential™ Silver & Gold packages
 - When challenges arise – the most efficient & cost-effective way



One stop service at GenScript: gene synthesis → Subcloning → PROtential™ → Scale up protein production

GenScript's Solution for Expression Optimization



Name	Service Type	Expression System(s)	Price	Timeline	Description
Standard	Protein Expression Evaluation	<i>E. coli</i> (SC1653-B)	Starting from \$280	1-2 weeks	Test 1 condition for soluble expression in a customer chosen bacterial strain
		Insect (SC1653-I)	Starting from \$400	3-4 weeks	Test 1 condition for soluble expression in Sf9, Sf21, S2 or Hi-5 cells
		Mammalian (SC1653-M)	Starting from \$500	3-4 weeks	Test 1 condition for soluble expression in CHO or 293 cells
		All 3 Systems (SC1653-3S)	Starting from \$950	3-4 weeks	Test soluble expression 1 condition/expression system (<i>E. coli</i> , Insect, Mammalian); total 3 systems
Silver	Protein Expression Optimization	<i>E. coli</i> (SC1667)	Login to inquire	2-3 weeks	<ul style="list-style-type: none"> • Test 8 different conditions • Optimize growth temperature, media components & inducer concentrations • Identify the best expression condition with your chosen vector and bacterial strain
Gold	Protein Expression Optimization	<i>E. coli</i> (SC1668)	Login to inquire	4-8 weeks	<ul style="list-style-type: none"> • Test 48 different conditions • Optimize growth temperatures, media components, inducer concentrations, promoters, host cell strains & fusion partners • Robust, industry-first, high-throughput expression and solubility optimization matrix

Case Study- Protein Expression Optimization



◆ Challenges:

- Pilot purification (final yield 1mg/L)- ~ 28kDa protein;
- The protein can be only purified from the soluble part
- 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

Case Study- Protein Expression Optimization



Strategy execution: HT expression testing

	1	2	3	4	5	6
A	X-1	X-2	X-3	X-4	X-5	X-6
B	X-7	X-8	X-9	X-10	X-11	X-12
C	X-13	X-14	X-15	X-16	X-17	X-18
D	X-19	X-20	X-21	X-22	X-23	X-24
E	X-25	X-26	X-27	X-28	X-29	X-30
F	X-31	X-32	X-33	X-34	X-35	X-36
G	X-37	X-38	X-39	X-40	X-41	X-42
H	X-43	X-44	X-45	X-46	X-47	X-48

Conditions X-1 to X-48

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 Choose Expression System
- 2 Optimize Protein Expression
- 3 Protein Refolding
- 4 Protein Purification
- 5 Difficult-to-Express Proteins
- 6 Advantages of GenScript Protein Services

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.

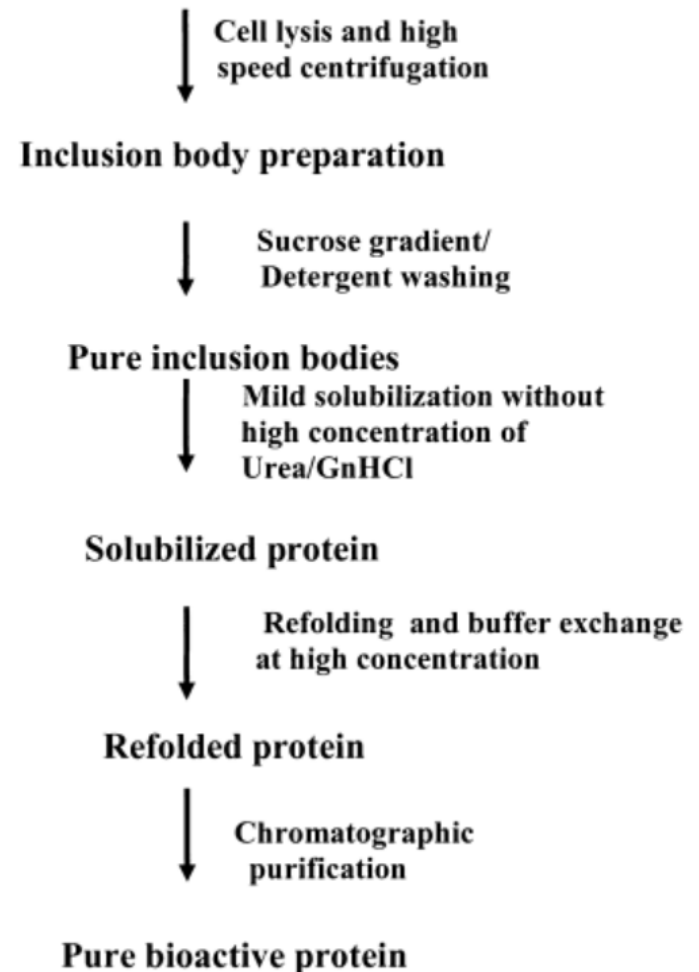
◆ Benefit

- 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

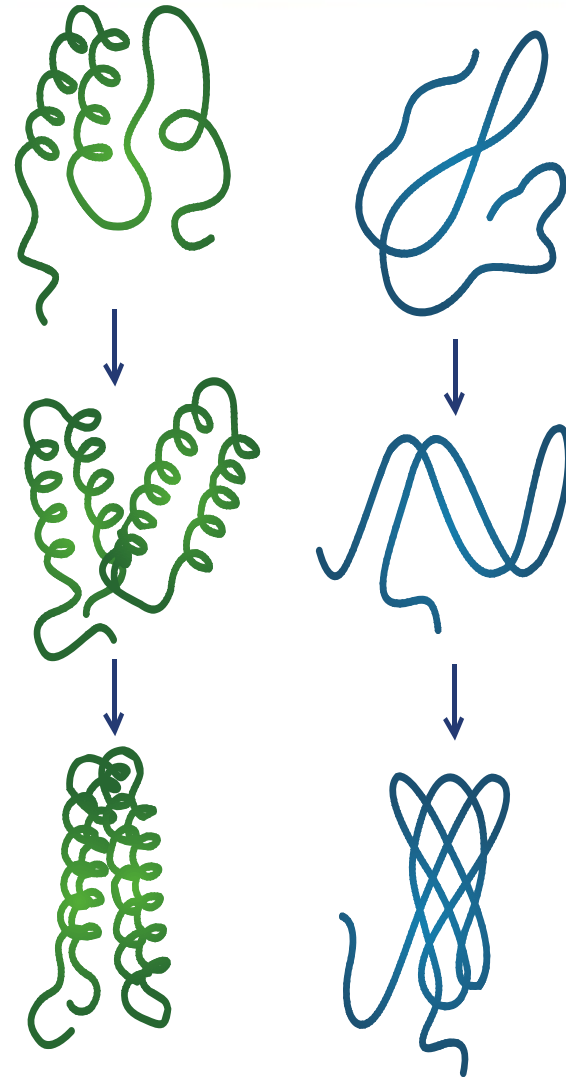
E. coli expressing protein as inclusion bodies



Protein Refolding Introduction



- ◆ All the information necessary for folding the peptide chain into its native structure is contained in the primary amino acid sequence of the peptide.
- ◆ The native form of a protein has the thermodynamically most stable structure.
- ◆ 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.



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.



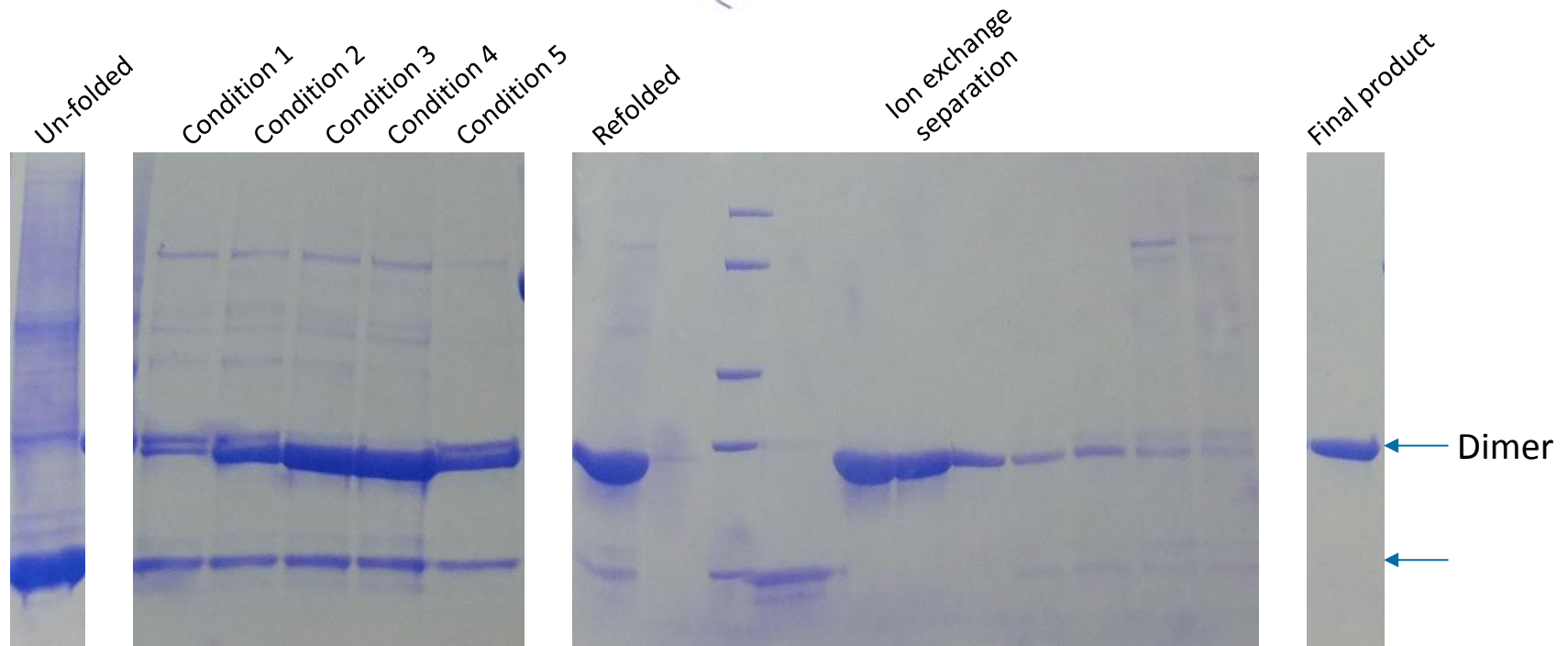
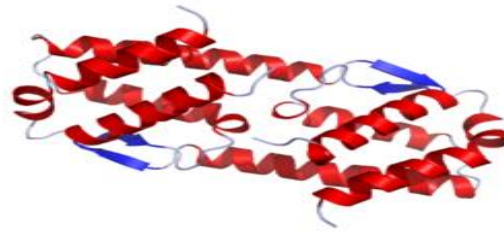
- ◆ Refolding conditions must be optimized for each individual protein.
- ◆ Important variables are:
 - buffer type
 - pH
 - ionic strength
 - Additives, often in combination (glycerol, redox reagents, saccharides, amino acids, metal ion, detergents, chaperones)

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

Case Study: Protein Refolding



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





- 1 Choose Expression System
- 2 Optimize Protein Expression
- 3 Protein Refolding
- 4 Protein Purification
- 5 Difficult-to-Express Proteins
- 6 Advantages of GenScript Protein Services

Protein Purification



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



Affinity Chromatography



◆ Affinity tags

- His
- GST
- MBP
- Flag
- SUMO

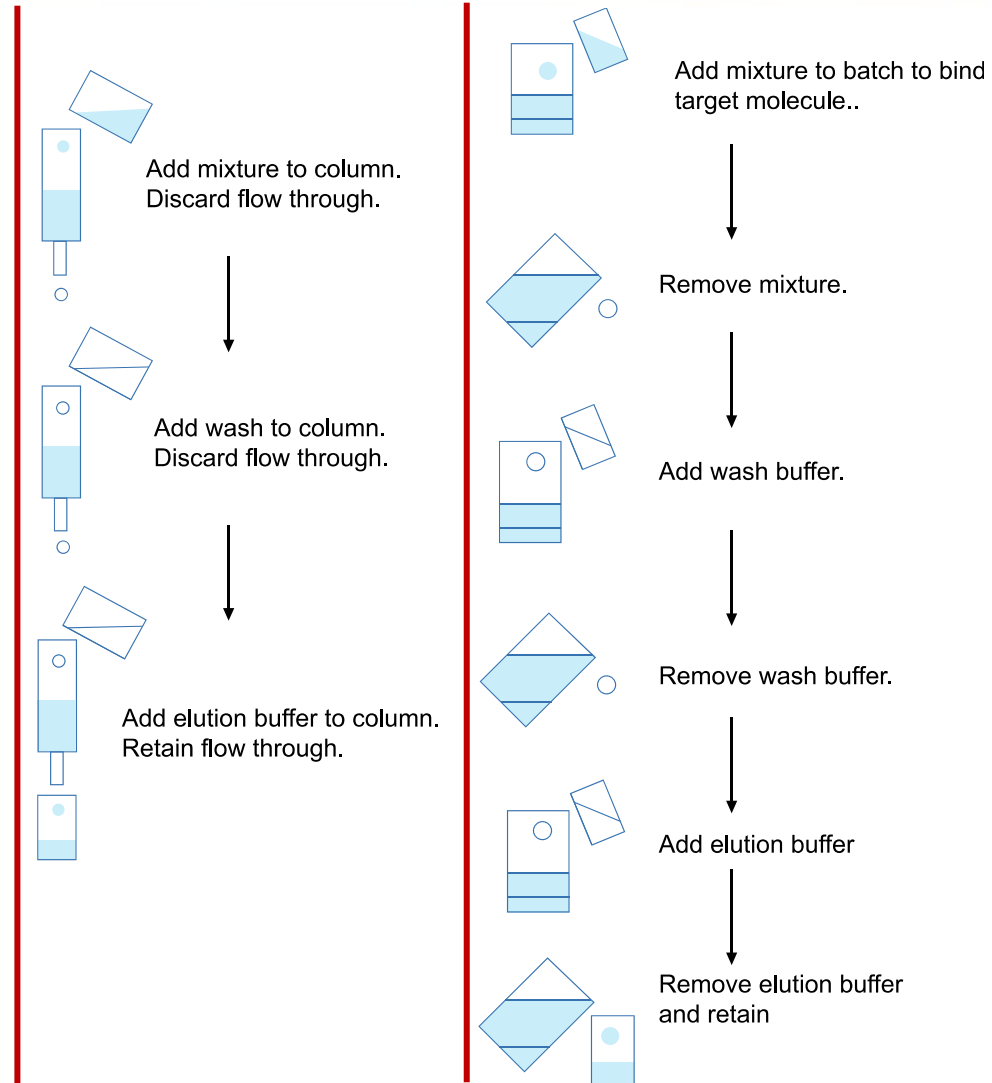
◆ Matrices & elution conditions

◆ Methods

- Column
- Batch

◆ Tag removal by site-specific protease

- TEV
- Enterokinase
- SUMO protease



Endotoxin Removal



◆ What are endotoxins?

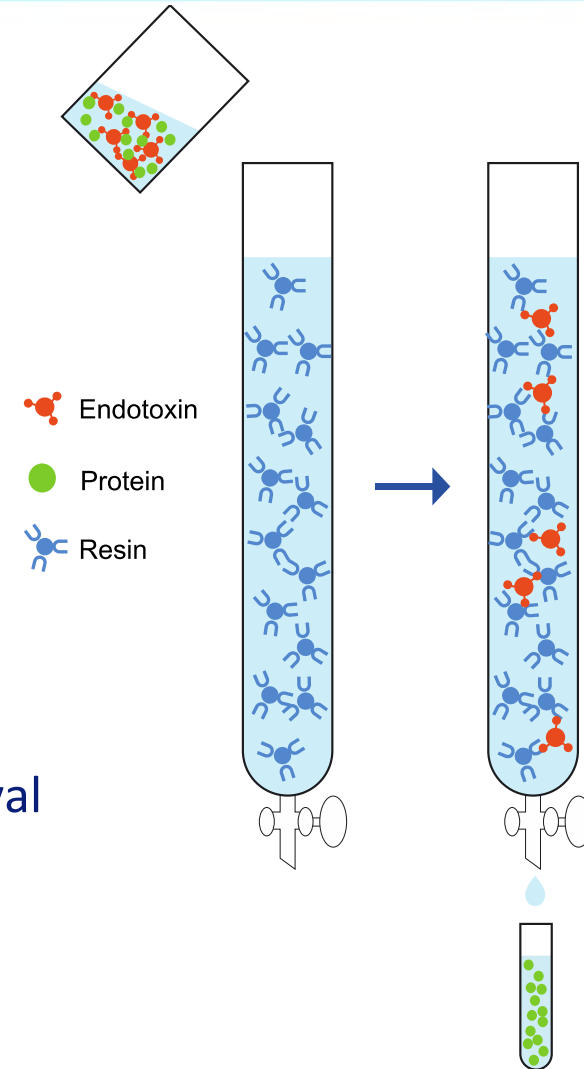
- Endotoxins, also known as lipopolysaccharides (LPS), are large molecules found in the outer membrane of Gram-negative bacteria, which elicit strong immune responses in animals.

◆ 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

GenScript Endotoxin Removal Services: Endotoxin ≤ 1 EU/ μ g; ≤ 0.1 EU/ μ g; ≤ 0.01 EU/ μ 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/ μ g	0.02 – 0.04 EU/ μ g	0.016 – 0.032 EU/ μ g

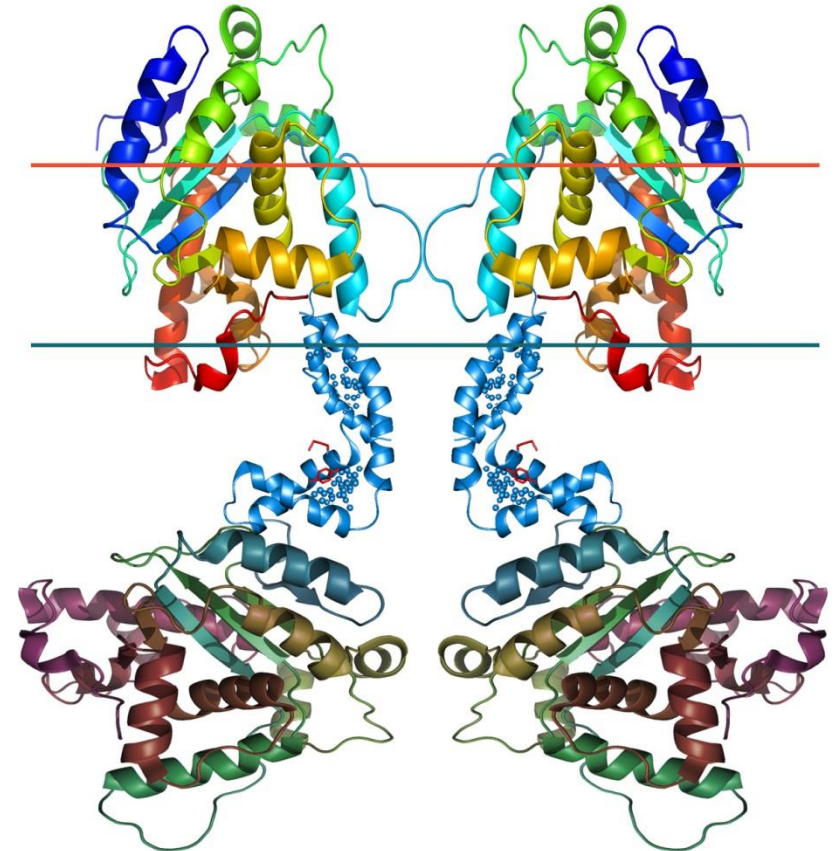


- 1 Choose Expression System
- 2 Optimize Protein Expression
- 3 Protein Refolding
- 4 Protein Purification
- 5 **Difficult-to-Express Proteins**
- 6 Advantages of GenScript Protein Services

Difficult-to-Express Proteins



- ◆ Membrane protein
- ◆ Toxic protein
- ◆ Prone-to-degrade protein



Membrane Protein



Definition: Membrane proteins are proteins that interact with biological membranes.

- Targets of over 50% of all modern medicinal drugs.
- 20-30% of all genes in most genomes encode membrane proteins.

❖ Expression

- Insect cells or mammalian cells;
- Budded baculovirus or virus-like particle;
- Cell-free

❖ Purification

- Detergent screen;
- Phosphate lipid ;
- Nano-disc;

Membrane Protein Expression Work Flow



Gene synthesis and subcloning into transfer vector

Generating the Recombinant Bacmid DNA

Transfecting Insect cells

Amplifying Baculovirus

Expression optimization

Large scale expression and harvest cell pastes

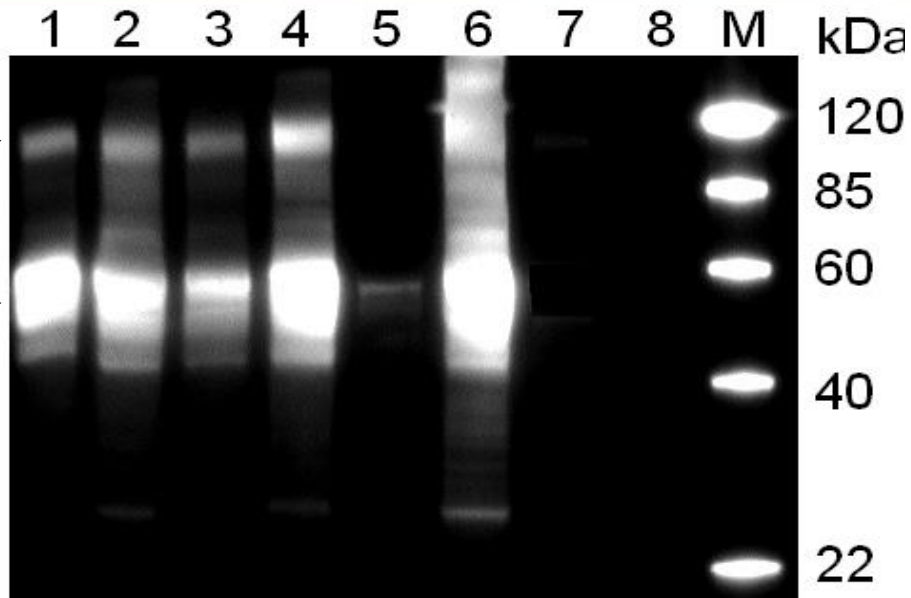
Homogenizing cells with a glass-Teflon tissue homogenizer

Centrifugation at low speed

Centrifugation at ultra speed and re-suspend the resulting pellet

- ◆ Deliver recombinant Baculovirus, or cell pastes, or prepared membrane protein to the customer.
- ◆ Deliver functional budded virus containing membrane protein to the customer.
- ◆ Provide high throughput screening of membrane proteins and deliver expression evaluation report to the customer.

Case Study: Membrane Protein



- ❖ 25 L expression of GPCR,
Expression Level: 30-40 mg/L
- ❖ Deliver the cell pastes in 6 weeks.

WB analysis of the membrane preparation

1. Whole cell lysate of GPCR before sonication
2. Whole cell lysate of GPCR after sonication
3. Supernatant after centrifugation of cell lysate for 10 minutes at 8k rpm
4. Pellet after centrifugation of cell lysate for 10 minutes at 8k rpm
5. Supernatant after ultracentrifugation of cell lysate for 45 min at 42k rpm
6. Pellet after ultracentrifugation of cell lysate for 45 min at 42k rpm
7. Negative control: supernatant after centrifugation of Sf9 cell lysate for 10 minutes at 8k rpm
8. Negative control: pellet after centrifugation of Sf9 cell lysate for 10 minutes at 8k rpm

Antibody: anti-His monoclonal-antibody (Genscript, Cat.No. A00186)



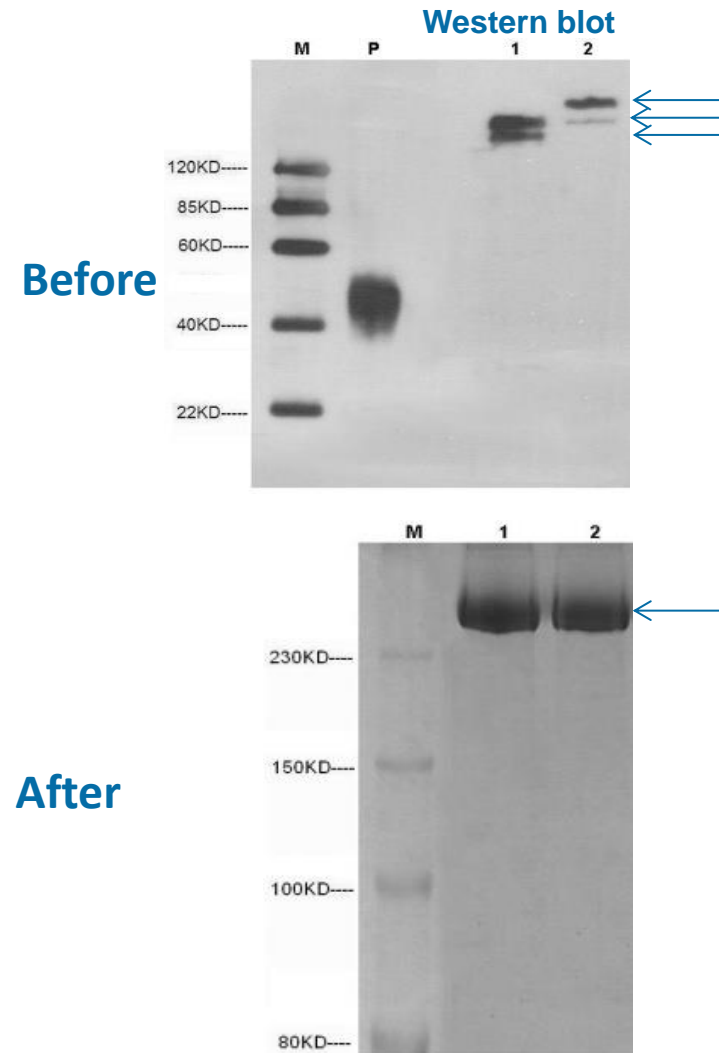
- ◆ Definition: Toxic proteins defined here as proteins that cause cell death or severe cultivation and maintenance defects during the growth phase when their genes were introduced into E. coli strain.
 - Mostly due to leaking expression
 - ~80% protein growth and expression problems are caused by the toxicity of proteins

- ❖ Strategies in solving the problem
 - Promoter selection
 - Suppress basal expression from leaky inducible promoters
 - Tight control of plasmid copy numbers
 - Protein production as inactive (insoluble) forms

Case Study: Prone-to-Degrade Protein



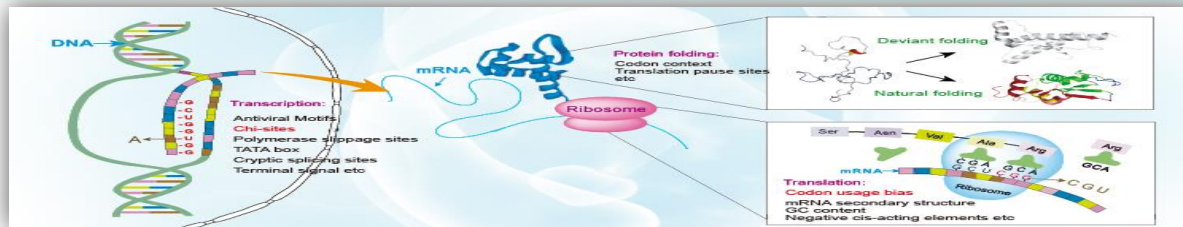
- ◆ Inconsistency of measured concentration
- ◆ Trouble shooting:
 - Transfection methods
 - Cell lysis
 - Purification
- ◆ Challenge: DNase is only partially responsible for the protein degradation. This protein itself is prone-to-degrade.
- ◆ Solutions:
 - Remove DNase
 - Add protease inhibitor to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





- 1 Choose Expression System
- 2 Optimize Protein Expression
- 3 Protein Refolding
- 4 Protein Purification
- 5 Difficult-to-Express Proteins
- 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 & Stable cell lines

Large Scale Protein Production

Upstream & Downstream Process
development, fermentation, GLP-compliant
Bioprocess Services

High Throughput Protein Variants

Largest high-throughput capacity in the industry,
proprietary platforms, 30 days for 1,000 protein
variants

Other Protein Services

Endotoxin removal, codon optimization, custom
purification, protein characterization, refolding

Protein Expression Evaluation & Optimization

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

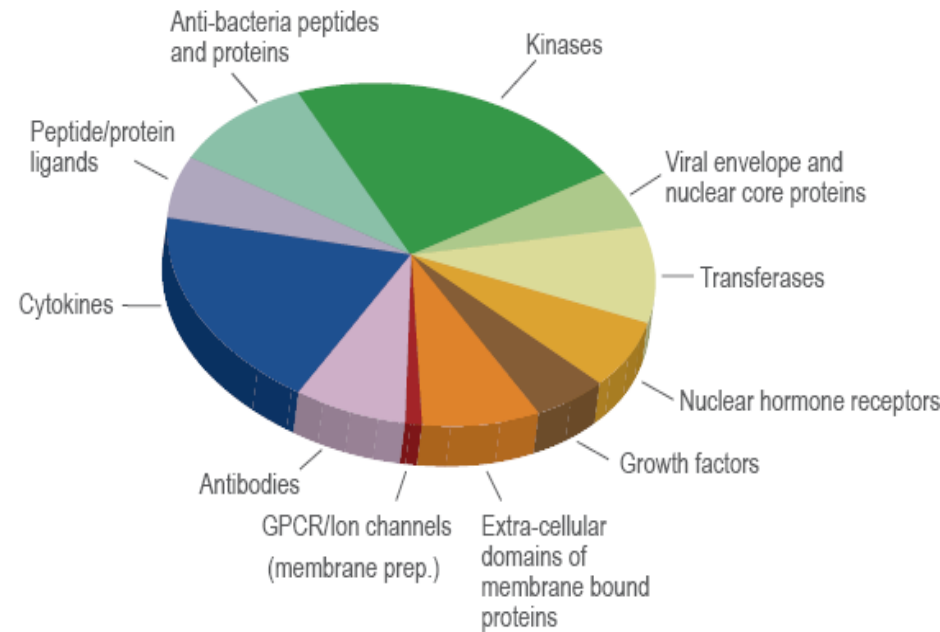
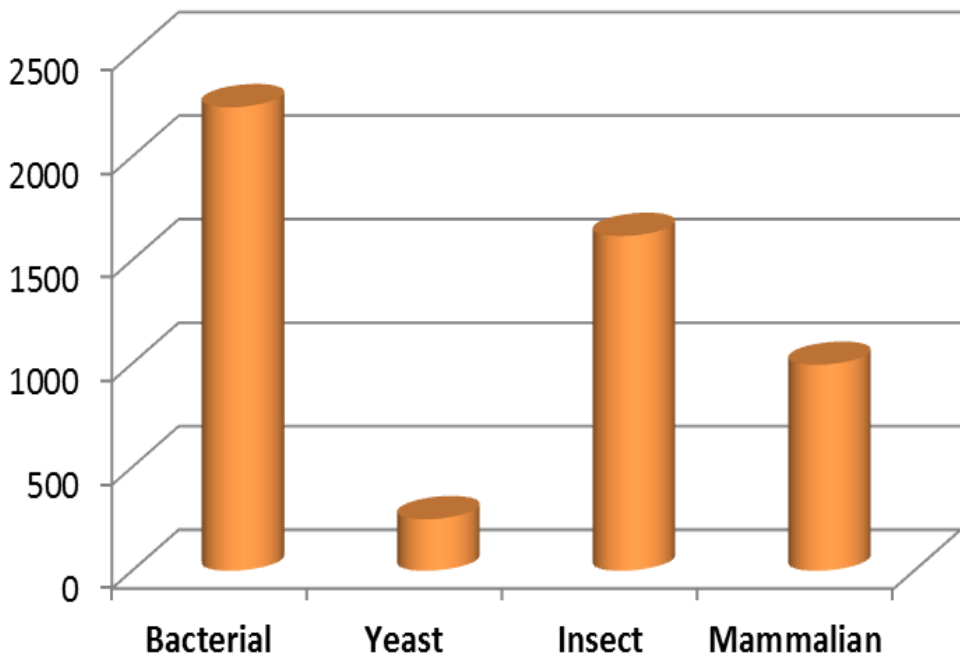
Chemical Protein Synthesis

Alternative method to produce high purity
functional proteins for hard-to-express proteins

Structural Biology

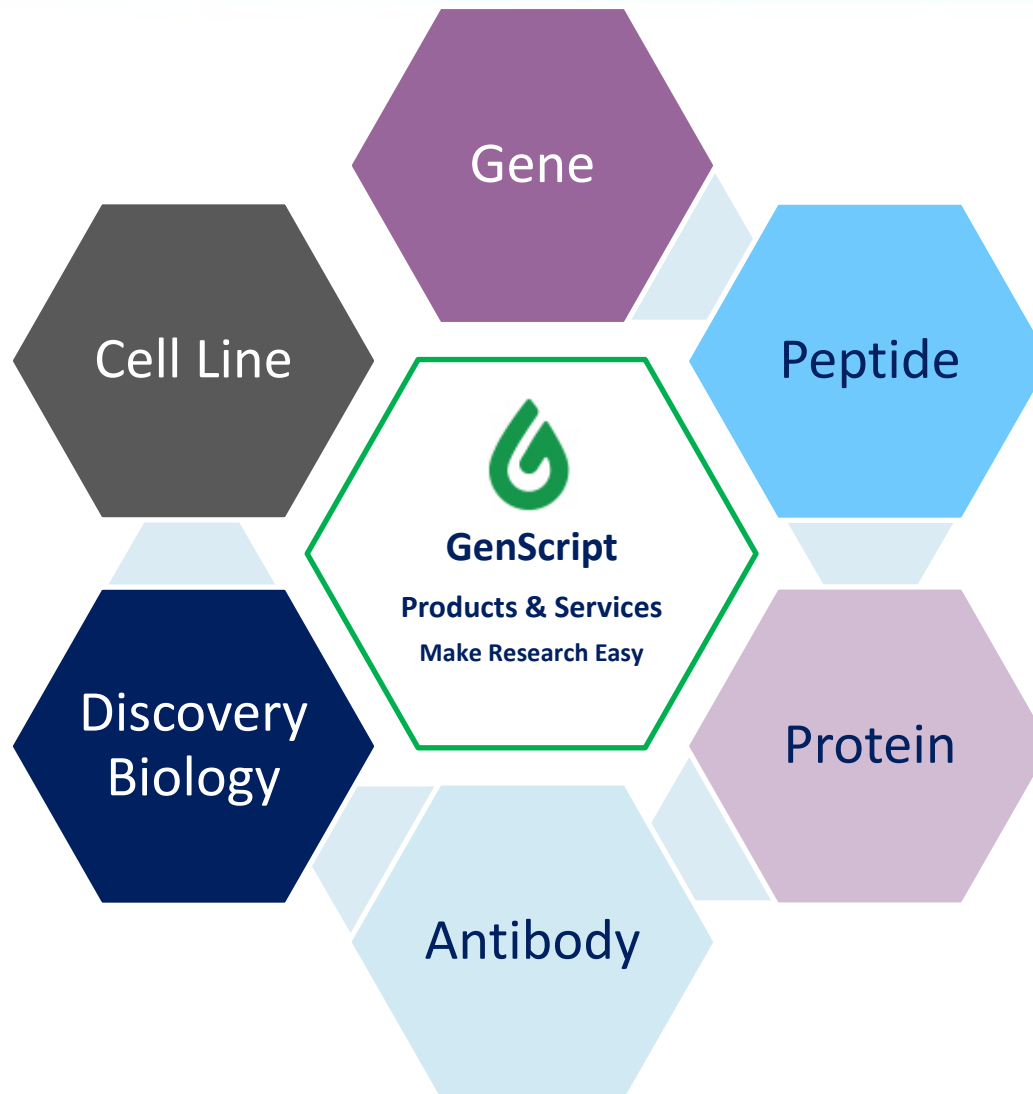
CrystalPro™ Gene-to-Structure Services, high
purity protein preparation, crystal, co-crystal
structure determination

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>



June 18, 2014/ 2:00 pm EST

Can CRISPR/Cas9 off-target genomic editing be avoided? Ways to improve target specificity - *Maxine Chen, Ph.D*



June 25, 2014/ 2:00 pm EST

Building a Synthetic Eukaryotic Genome – Sc2.0 - *Leslie Mitchell, Ph.D., NYU Langone Medical Center*

Promo Code: **BacPowerUP**

- Free additional amount of protein from 3mg to 5mg for any BacPower™ Guaranteed Packages
- Valid till 8/15/2014