

Codon optimization: Why & how to design DNA sequences for optimal soluble protein expression

Rachel Speer, Ph.D.



Codon optimization: Why & how to design DNA sequences for optimal soluble protein expression



- 1 Why do codons matter?
- 2 OptimumGene principles & performance
- 3 Case Studies
- 4 How to get better protein expression

Protein Expression Overview



Select/Design the end product
(amino acid sequence)



Choose expression system



Design expression clone
(DNA construct)



Express the protein

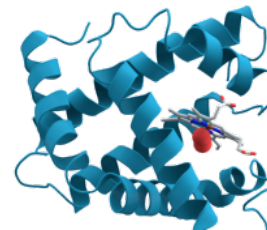
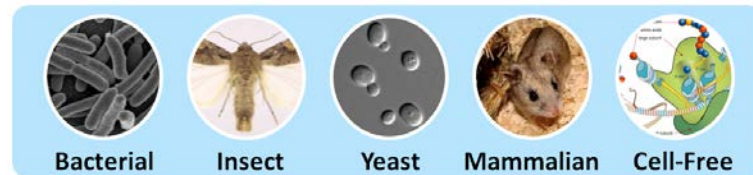


Purify the protein



Characterize the protein

MGVHECPAWLWLLLSLLSLPLGLPVLGAPPRLIC...



Protein Expression Overview



Select/Design the end product
(amino acid sequence)



Choose expression system



Design expression clone
(DNA construct)



Express the protein

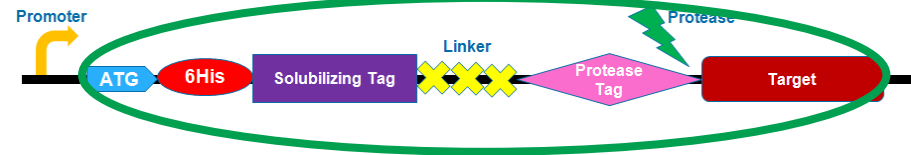


Purify the protein



Characterize the protein

Codon Optimization



Why do Codons Matter? The Facts



- ◆ Redundancy in the genetic code
- ◆ Synonymous mutations affect protein expression rates up to 1000-fold.
- ◆ Synonymous mutations can also alter protein conformation, PTM, stability, and function.

		Second Letter					
		U	C	A	G		
1st letter	U	UUU Phe UUC UUA Leu UUG	UCU Ser UCC UCA UCG	UAU Tyr UAC UAA Stop UAG Stop	UGU Cys UGC UGA Stop UGG Trp	U C A G	3rd letter
	C	CUU CUC Leu CUA CUG	CCU CCC Pro CCA CCG	CAU His CAC CAA Gln CAG	CGU CGC Arg CGA CGG	U C A G	
	A	AUU AUC Ile AUA AUG Met	ACU ACC Thr ACA ACG	AAU Asn AAC AAA Lys AAG	AGU Ser AGC AGA Arg AGG	U C A G	
	G	GUU GUC Val GUA GUG	GCU GCC Ala GCA GCG	GAU Asp GAC GAA Glu GAG	GGU GGC Gly GGA GGG	U C A G	

Codon Optimization:

Introducing synonymous mutations that favor efficient soluble protein expression

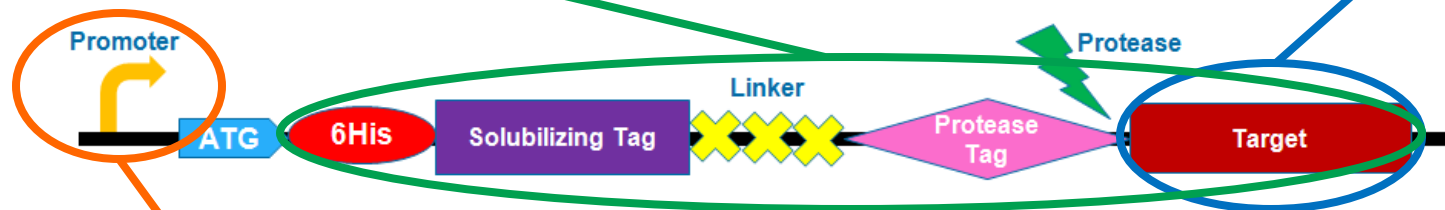
Optimized	AGTTTTCAGGTGAGGTCCGCCCGTT
Original	AGCTTCCCGGGATGAGGGCCCCCGGTT

What Codon Optimization is – and isn't



Codon Optimization:
Introducing synonymous mutations that favor efficient soluble protein expression

Protein Design:
changing the amino acid sequence



Expression strategy:
Selecting promoter, tags,
etc.



Observations:

- Species-specific bias in codon use and tRNA abundance
- Heterologous protein expression is often inefficient

Theory:

- Rare codons reduce protein expression

Solutions:

- Express tRNA to remove bias in the host cells
- Alter the gene to replace rare codons with preferred ones:
 - site-directed mutagenesis
 - *de novo* gene synthesis with codon optimization

Codon Optimization Significantly Increases Protein Expression Levels



Codon optimization can improve expression of human genes in *Escherichia coli*: A multi-gene study.

Burgess-Brown NA *et al. Protein Expr Purif.* May 2008; 59(1): 94-102

Gene Name	Native			Synthetic			Expression	Solubility
	1	2	3	1	2	3	Syn vs Nat	Syn vs Nat
CBR1	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	▲	▲
CBR3	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	▲	▲
GMD5	Not Expressed	Not Expressed	Not Expressed	Not Expressed	Expressed, Soluble and Purified	Expressed, Soluble and Purified	▲	▲
HADH2	Not Expressed	Expressed	Expressed	Not Expressed	Expressed, Soluble and Purified	Expressed	▲	▲
HSD17B2	Expressed	Expressed	Not Expressed	Expressed	Expressed	Not Expressed	▲	
HSD17B4	Not Expressed	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	▲	▲
MGC4172	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified		
PECR	Not Expressed	Not Expressed	Not Expressed	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed	▲	▲
RETSR2	Expressed	Expressed, Soluble and Purified	Expressed	Expressed	Expressed	Expressed	▲	
SPR	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified	Expressed, Soluble and Purified		

Expressed
Expressed, Soluble and Purified
Not Expressed

▲ Targets shown improvement of expression and/or solubility with synthetic gene after codon optimization

1. Total Cellular Protein
2. Soluble Fraction
3. Eluted Fraction

Codon Adaptation is not the most important factor for protein yield



Coding-sequence determinants of gene expression in *Escherichia coli*.

Kudla G, Murray AW, Tollervey D, Plotkin JB. *Science*. 2009 Apr 10;324(5924):255-8.

- 154 synthetic GFP genes with random synonymous mutations
- 250-fold variation in fluorescence
- 44% of variation explained by 5' mRNA free energy (nt -4 to +37)

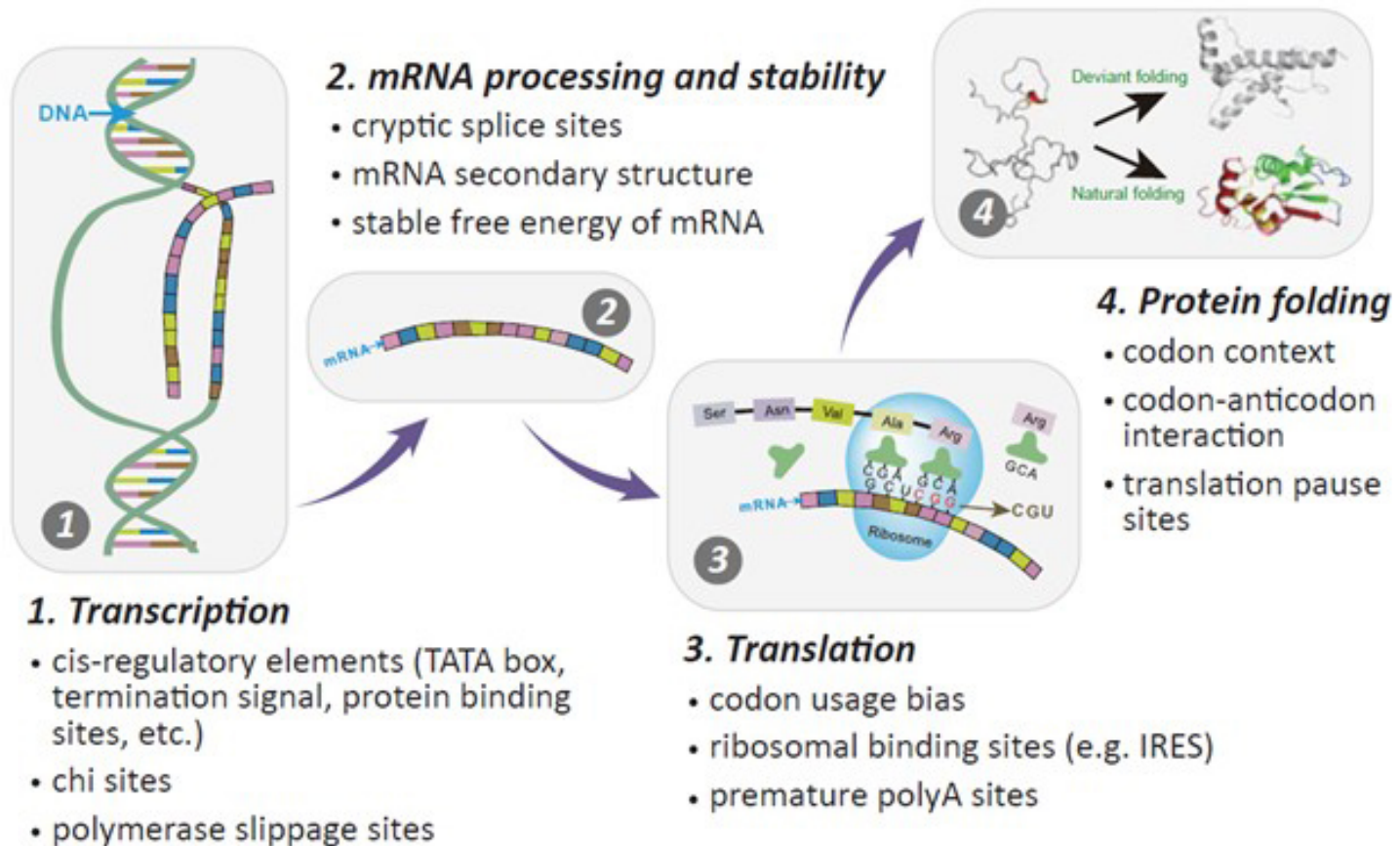


The anti-Shine-Dalgarno sequence drives translational pausing and codon choice in bacteria.

Li GW, Oh E, Weissman JS. *Nature*. 2012 Mar 28;484(7395):538-41.

- Variation in Translation Rate does not correlate with rare codon use
- Orthogonal ribosomes with altered anti-SD sequences: pausing results from hybridization between 16s rRNA and SD-like sequences in mRNA

Many sequence features influence protein expression



Evidence-Based Codon Optimization: OptimumGene



Transcriptional Efficacy:

- GC content
- CpG dinucleotides content
- Cryptic splicing sites
- Negative CpG islands
- SD sequence
- TATA boxes
- Terminal signal

Translation Efficiency:

- Codon usage bias
- GC content
- mRNA secondary structure
- Premature PolyA sites
- RNA instability motif (ARE)
- Stable free energy of mRNA
- Internal chi sites and ribosomal binding sites

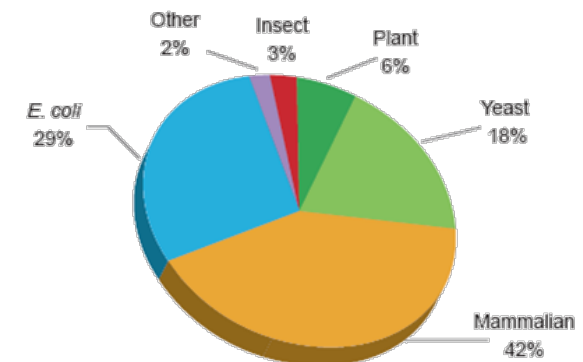
Protein Refolding:

- Codon usage bias
- Interaction of codon and anti-codon
- Codon-context
- RNA secondary structures

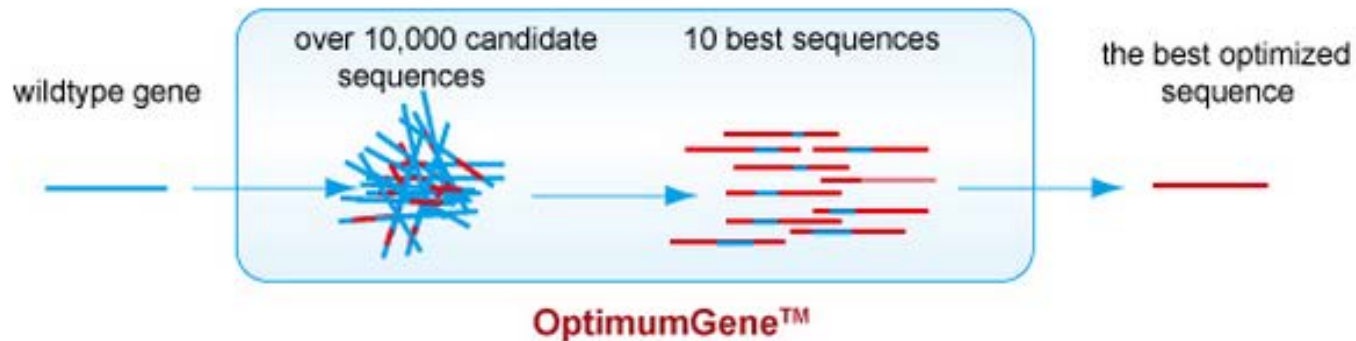
Flexibility to adjust the weight of different factors or add customized constraints:

- Filter out restriction sites
- Reduce similarity between library members
- Alternative codon tables / condition-specific codon preferences

GenScript has optimized over **50,000 sequences** in all major expression systems.



Patented Bioinformatic Algorithm powers OptimumGene



Liu *et al.* **Method of sequence optimization for improved recombinant protein expression using a particle swarm optimization algorithm.** US Patent 8,326,547, issued December 4, 2012.

OptimumGene™ Improves Protein Expression Better than Competitors' Optimization

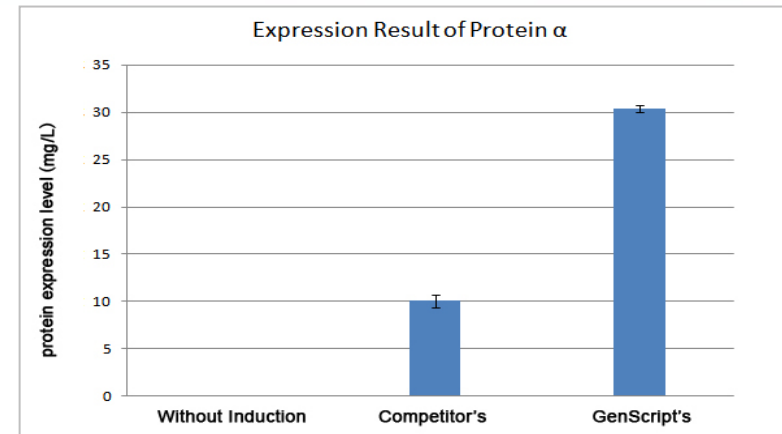
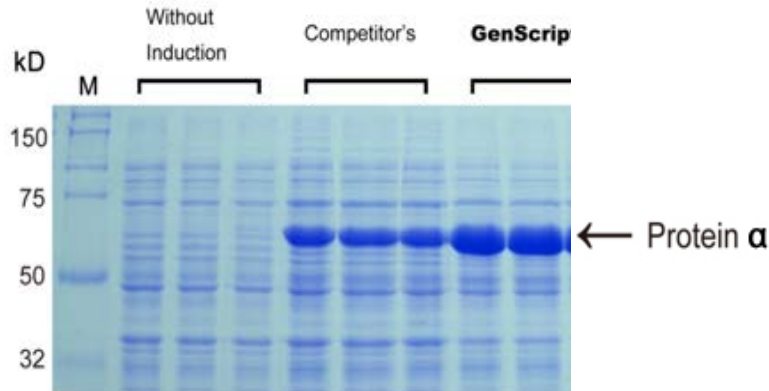


Fig. 1: Expression Result of Protein α after Codon Optimization. The expression level of Protein α using GenScript's OptimumGene™ Codon Optimization is **3** times more than that of competitor's.

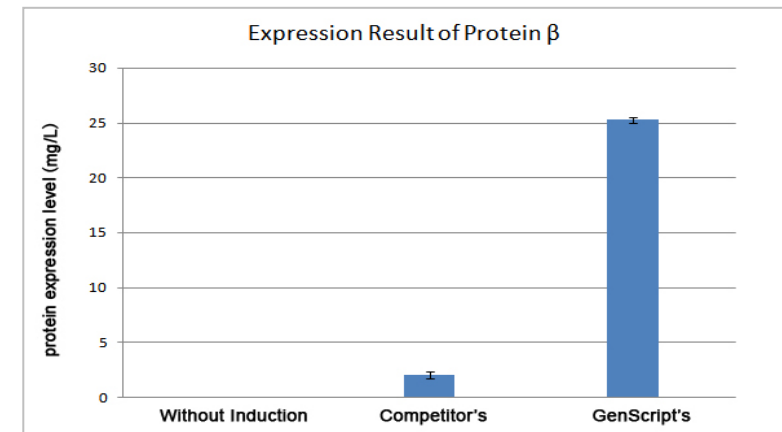
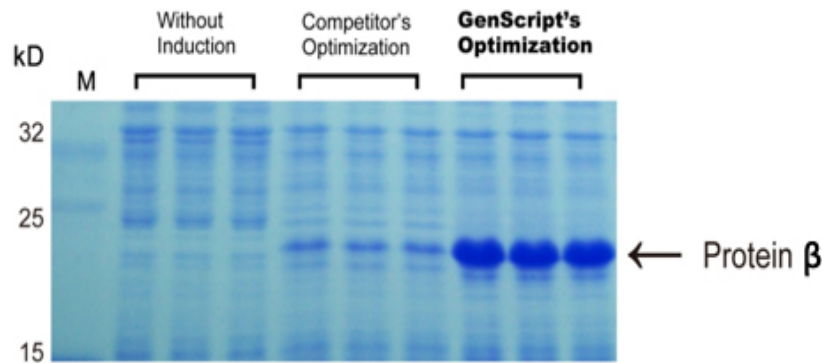
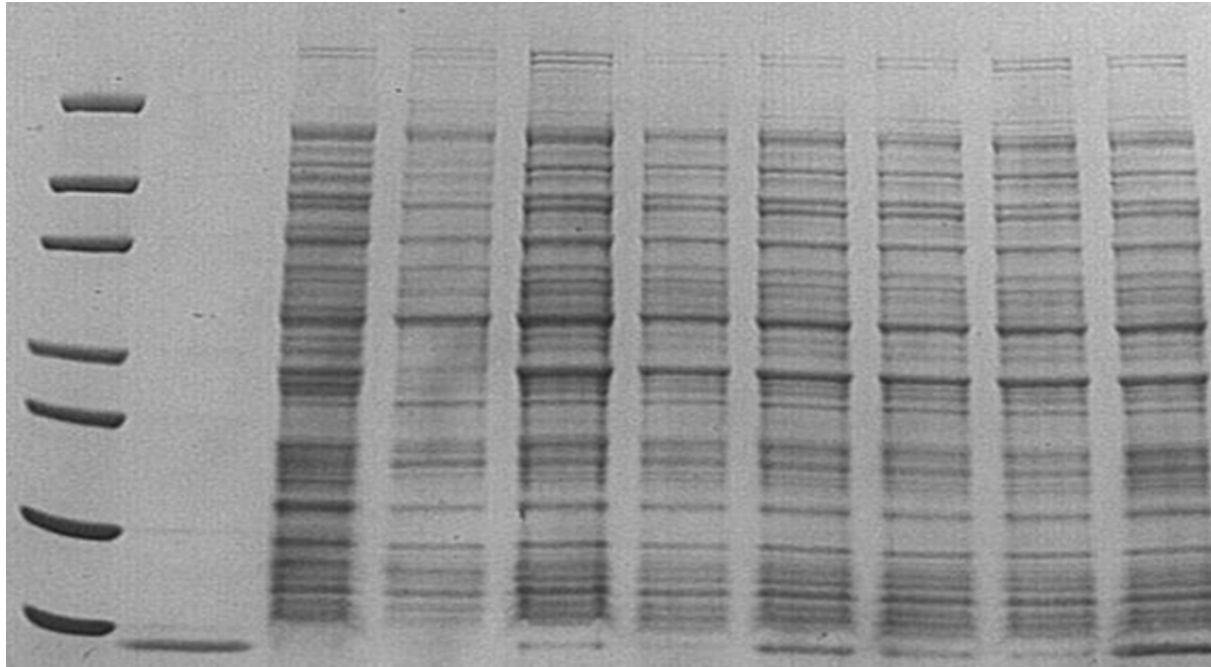


Fig. 2: Expression Result of Protein β after Codon Optimization. The expression level of Protein β using GenScript's OptimumGene™ Codon Optimization is **13** times more than that of competitor's.

We routinely test and improve our algorithm



Lane 1 2 3 4 5 6 7 8 9 10



Lane	1	2	3	4	5	6	7	8	9	10
Sequence	MW marker	Purified hIGF-1 (PC)	BL21 cell lysate (NC)	WT (non-optimized) hIGF-1	Opti-0	Opti-1	Opti-2	Opti-3	Opti-4	Opti-5
Yield (mg/L)	--	--	--	Not detectable	7.7	3.1	18.5	11.4	5.4	28.5

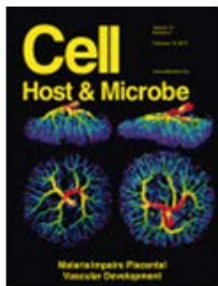
Hundreds of papers cite GenScript for codon-optimized gene synthesis



“Humanization and optimization of codon usage was performed (GenScript) owing to **poor expression of the original zebrafish lyn in HEK293 cells.**”



“...IFP1.4 gene was de novo synthesized by **GenScript Company**, based on the available protein sequence. The DNA sequence was **optimized with proprietary OptimumGene algorithm (GenScript)**...”



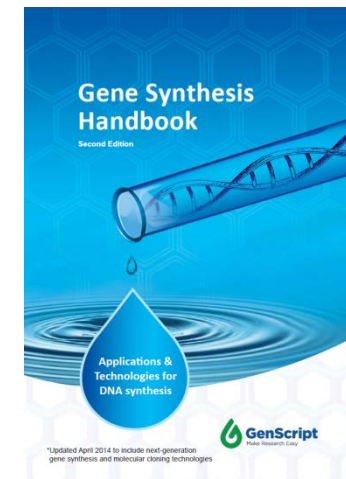
“...The following genes were **codon optimized and synthesized (Genscript)**:...”

[Resources](#) » [Reference Databases](#) » [Citations Database](#)

By Category

- ▼ Peptide Services(1862)
- ▼ Antibody Engineering(1)
- ▼ Animal Model Services(4)
- ▼ Bio-Assay Center(3)
- ▼ Cell Line Services(6)
- ▼ Antibody Services(628)
- ▼ Gene Services(3169)
- ▼ Protein Services(160)
- ▼ Bioinformatics Tools(332)
- ▼ Catalog Products(3042)

[View all \(10456\)](#)



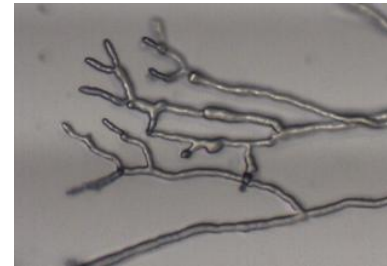
Case Study 1: Neurospora Circadian Clock



nature

Non-optimal codon usage affects expression, structure and function of clock protein FRQ

Zhou M. *et al. Nature*. 2013 Mar;495 (7439); 111 – 5



Codon Optimization performed in only specific regions of the gene:

“...Optimized frq sequences (synthesized by Genscript) ...In the m1-frq construct, only the codons upstream of the predicted intron branch point were optimized as m-frq. For the m2-frq construct, only the codons downstream of the intron 3' end were optimized as m-frq...”

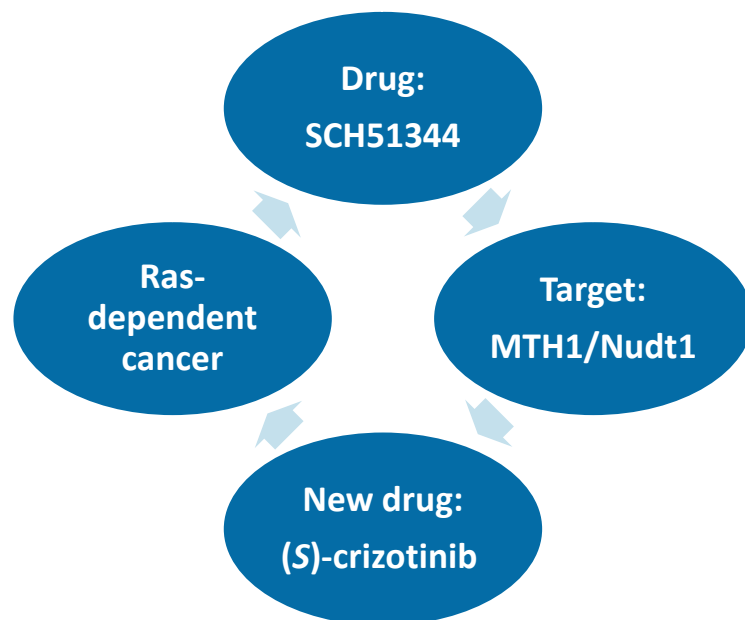
Case Study 2: Crystal Structure & PK/PD for Cancer Drug Discovery



nature

Stereospecific targeting of MTH1 by (S)-crizotinib as anticancer strategy

Huber KV, et al. *Nature*. 2014 Apr 10;508(7495):222-7.



- Codon-optimized gene synthesis from GenScript was used to express Nudt1 for enzymatic assays and crystallization studies.

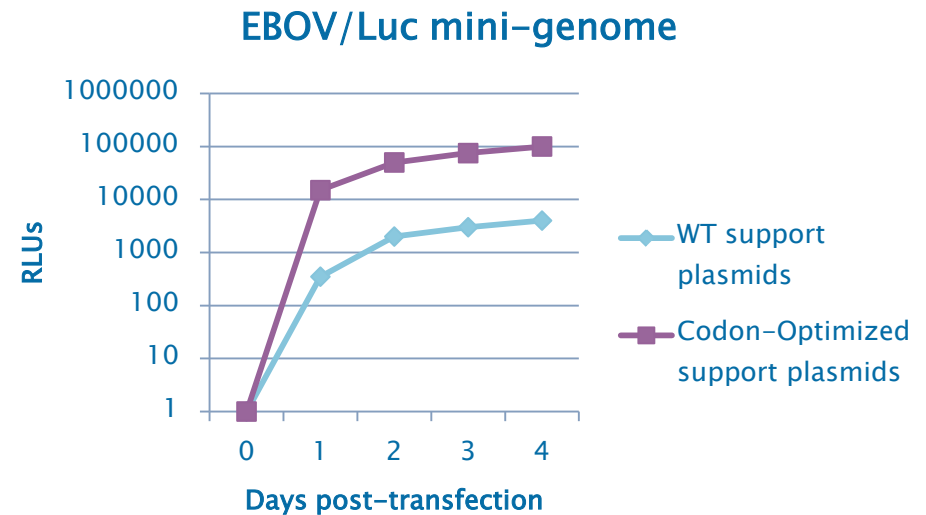
Case Study 3: Antiviral Drug Discovery for Ebola Virus



High-throughput, luciferase-based reverse genetics systems for identifying inhibitors of Marburg and Ebola viruses.

Uebelhoer *et al. Antiviral Res.* 2014 Jun;106:86-94.

- Codon-optimized EBOV gene provided by GenScript
- Codon-optimized support plasmids increased signal 2000-fold



Case Study 4: Combinatorial Library for Human Genome Editing



nature
biotechnology

A library of TAL effector nucleases spanning the human genome.

Kim Y, et al. *Nat. Biotechnol.* **31**, 251–258 (2013).

Table I. TALE specificity code

RVD	Nucleotide
NI	adenine
HD	cytosine
NG	thymine
NN	guanine

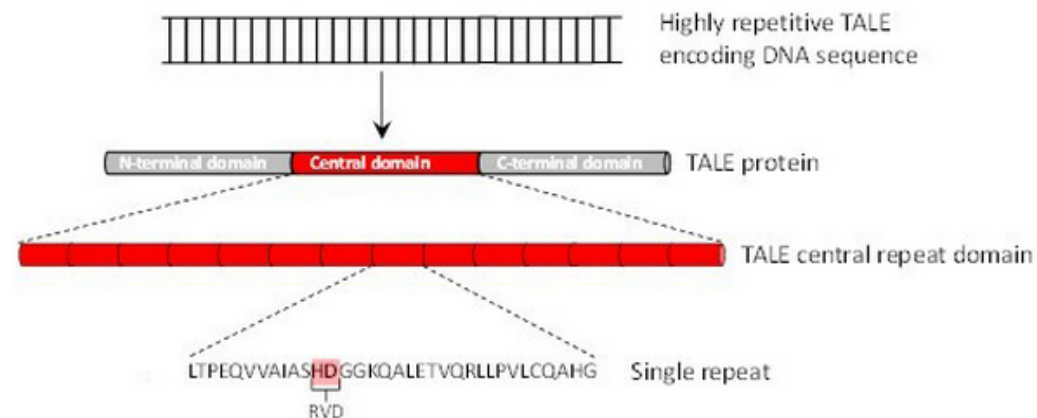
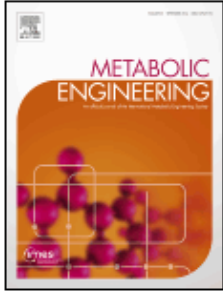


Fig 2. TALE protein organization

•Codon Optimized Gene Synthesis from GenScript was used to

1. Limit sequence similarity
2. Exclude rare codons
3. Guarantee accuracy of highly-repeated sequences

Case Study 5: Metabolic Engineering



Substantial improvements in methyl ketone production in *E. coli* and insights on the pathway from in vitro studies.

Goh EB *et al. Metab Eng.* 2014 Sep 18;26C:67-76.



- **Codon Optimized Gene Synthesis from GenScript** was used to
 1. improve metabolic pathway efficiency (\uparrow substrate influx, \downarrow diversion)
 2. improve GC content of gene from *M. luteus*, whose genome is 73% GC

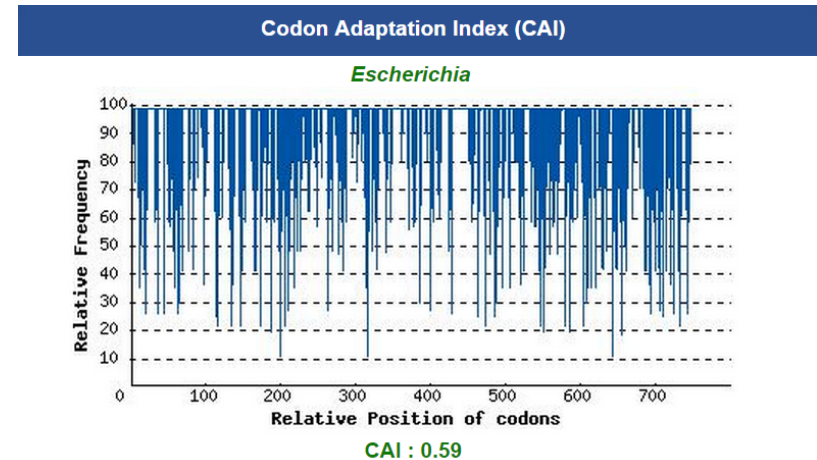
How to get codon-optimized genes



- Online Tools to identify rare codons



- Rare Codon Analysis Tool
- Codon Frequency Tables





- Request Free Codon Optimization using OptimumGene

Request Free Codon Optimization



•Quick & easy online form

	Email: gene@genscript.com
	Phone: 1-877-436-7274 (Toll-Free)

Information

Verify

Quote

Order

VeriSign

Design your Gene

1. Gene Sequence Information

* Gene name (for your records only):

☐ 5' sequence: ?

* Sequence: ☒ DNA ☐ Protein

☐ 3' sequence:

[Batch Orders \(xls\)](#)

2. Codon Optimization

☒ Yes ☐ No

3. Gene Synthesis

☒ Standard Service ☐ Rush Service ☐ Fast Service

4. Cloning ?

☒ Standard Cloning ☐ Express Cloning ☐ Custom Cloning

Vector:

Other Requirements:

5. Plasmid Prep ?

gene name

Codon Optimization

(* Required fields)

*Host expression organism:

Secondary expression organism:

(Optional) Note: If the secondary expression host is not closely related to the primary host, the optimization may result in a sequence with compromised codon usage index for both hosts. Selecting only the primary host will generate the most optimum sequence for the specified expression organism.

Optimization start position: bp, end position: bp

ORF start position: bp, end position: bp

Restriction sites to avoid: (example: BamHI, HindIII(AAGCTT))

Restriction sites to keep: (example: BamHI, HindIII(AAGCTT))

Stop codon needed:

☒ Yes. Sequence:

☐ No

Review your Free Optimization Report



Codon Adaptation Index (CAI)

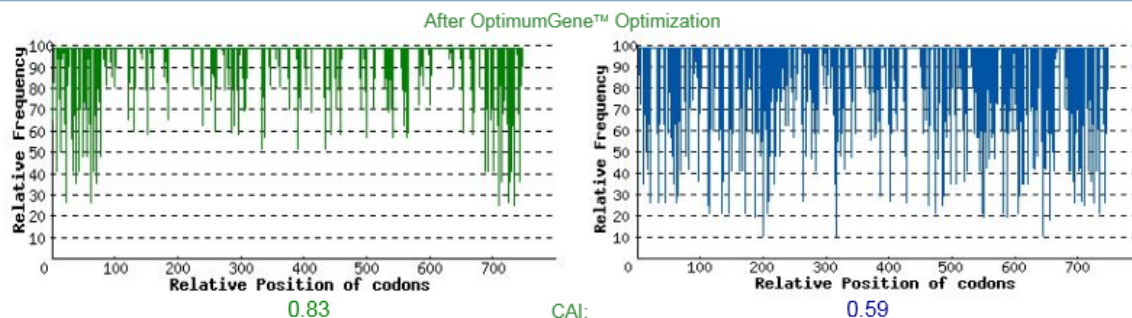


Figure 1a. The distribution of codon usage frequency along the length of the gene sequence. A CAI of 1.0 is considered to be perfect in the desired expression organism, and a CAI of > 0.8 is regarded as good, in terms of high gene expression level.

GC Content Adjustment

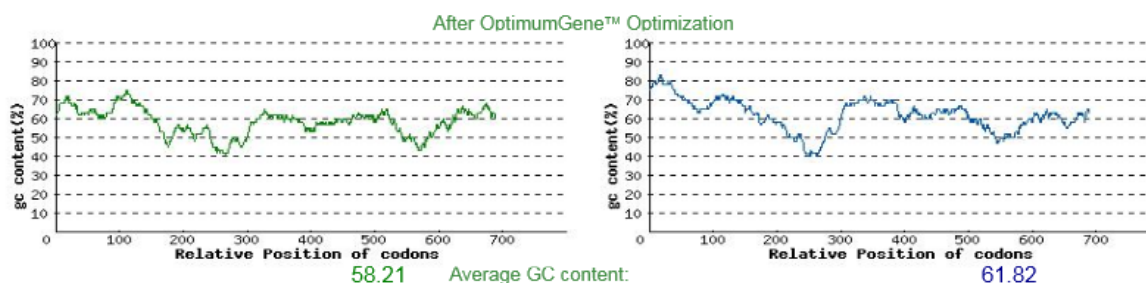
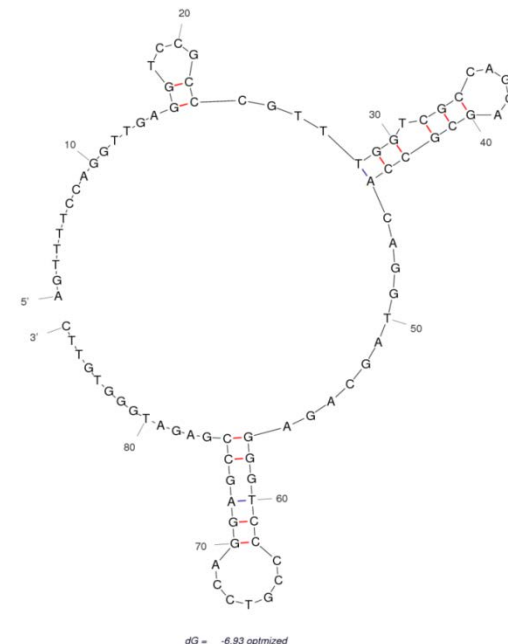


Figure 2. The ideal percentage range of GC content is between 30-70 %. Peaks of %GC content in a 60 bp window have been removed.



Helices in structure (all)

Helix	ΔG (kcal/mol)	Length	Position
1	-4.74	5	64-->68 ; 85<--81
2	-4.08	3	16-->18 ; 42<--40
3	-3.12	3	19-->21 ; 38<--36
4	-2.17	2	69-->70 ; 76<--75
5	-2.17	2	24-->25 ; 33<--32
6	-1.84	2	53-->54 ; 59<--58
7	-1.84	2	43-->44 ; 49<--48

Hairpins in structure (all)

Hairpin	ΔG (kcal/mol)	Length	Position
1	3.10	10	24-->...<--33
2	2.50	8	69-->...<--76
3	1.50	7	53-->...<--59
4	1.50	7	43-->...<--49

Order Gene Synthesis for your Codon-Optimized Gene



Recommended Services for your needs:	Low Price	Fast Turnaround	High-Volume	Long Genes
Custom Gene Synthesis Cat No. SC1010	✓ \$0.35/bp	✓ 9 business days	No min / max	≤8 kb
Rush Gene Synthesis Cat No. SC1575	Request a quote	✓ 4 business days	No min / max	≤2 kb
GenPlus™ High-Throughput Gene Synthesis Cat No. SC 1645	✓ \$0.23/bp	15 business days	✓ ≥25 genes	✓ ≤10 kb
GenPlus™ Economy Gene Synthesis Cat No. SC1681	✓ \$0.23/bp	25 business days	No min / max	✓ ≤10 kb
GenBrick™ Synthesis Cat No. SC1584	\$0.45/bp	23 business days	No min / max	✓ 8 - 15kb or more

GenScript Toolkit For Improving Protein Expression



**Select/Design the end product
(amino acid sequence)**



Choose expression system



**Design expression clone
(DNA construct)**



Express the protein



Purify the protein



Characterize the protein

GenPlus™ high-throughput gene synthesis
Gene Variant Library services

PROtential™ protein expression evaluation service

Codon-Optimized Gene Synthesis

BacPower™
YeastHIGH™
MamPower™
InsectPower™

FragPower™
Recombinant Antibody

FoldArt™ Refolding
ToxinEraser™ Endotoxin Removal

Protein Characterization Services

Related GenScript Webinars



Select/Design the end product
(amino acid sequence)



Choose expression system



Design expression clone
(DNA construct)



Express the protein



Purify the protein



Characterize the protein

Mutant library for protein engineering?
Combinatorial Library?
Truncation Variants?

Gene variant libraries: design, construction, and research applications



Presented by: Rachel Speer, Ph.D.
Originally aired May 21st and June 18th, 2014

On Demand

[View now](#)

Related GenScript Webinars



Select/Design the end product
(amino acid sequence)



Choose expression system



Design expression clone
(DNA construct)



Express the protein



Purify the protein



Characterize the protein

- Pros and Cons of different expression hosts
- Techniques for protein re-folding, protection from rapid degradation

Recombinant protein expression & purification: challenges and solutions



Presented by: Liyan Pang, Ph.D.
Originally aired June 11th and June 12th, 2014

On Demand

[View now](#)

Related GenScript Webinars



Select/Design the end product
(amino acid sequence)



Choose expression system



Design expression clone
(DNA construct)



Express the protein



Purify the protein



Characterize the protein

- Fusion partners/epitope tags
- e. coli strain, induction conditions, etc

Optimizing conditions for recombinant soluble protein production in *E. coli*



Presented by: Keshav Vasanthavada
Originally aired May 8th and June 24th, 2014

On Demand

[View now](#)

Related GenScript Webinars



Select/Design the end product
(amino acid sequence)



Choose expression system



Design expression clone
(DNA construct)



Express the protein



Purify the protein



Characterize the protein

- maximizing purity and yield

**Identify the optimal protein purification strategy
for your recombinant protein production**



Presented by: Liyan Pang, Ph.D.

November 12, 2014
8:00 am

[Register now](#)

November 12, 2014
2:00 pm

[Register now](#)

GenScript – The most cited biology CRO



References



- Belfield EJ, Hughes RK, Tsesmetzis N, Naldrett MJ, Casey R **The gateway pDEST17 expression vector encodes a -1 ribosomal frameshifting sequence.** *Nucleic Acids Res.* 2007;35(4):1322-32.
- Chu D *et al.* **Translation elongation can control translation initiation on eukaryotic mRNAs.** *EMBO J.* 2014 Jan 7;33(1):21-34.
- Goh EB *et al.* **Substantial improvements in methyl ketone production in E. coli and insights on the pathway from in vitro studies.** *Metab Eng.* 2014 Sep 18;26C:67-76.
- Huber KV, *et al.* **Stereospecific targeting of MTH1 by (S)-crizotinib as anticancer strategy.** *Nature.* 2014 Apr 10;508(7495):222-7
- Kim Y, *et al.* **A library of TAL effector nucleases spanning the human genome.** *Nat. Biotechnol.* **31**, 251–258 (2013).
- Kramer G, Boehringer D, Ban N, Bukau B. **The ribosome as a platform for co-translational processing, folding and targeting of newly synthesized proteins.** *Nat Struct Mol Biol.* 2009;16:589–597.
- Kudla G, Murray AW, Tollervey D, Plotkin JB. **Coding-sequence determinants of gene expression in Escherichia coli.** *Science.* 2009 Apr 10;324(5924):255-8. [Free Full Text](#)
- Li GW, Oh E, Weissman JS. **The anti-Shine-Dalgarno sequence drives translational pausing and codon choice in bacteria.** *Nature.* 2012 Mar 28;484(7395):538-41. [Free Full Text](#)
- Mellitzer A. *et al.* **Synergistic modular promoter and gene optimization to push cellulase secretion by Pichia pastoris beyond existing benchmarks.** *J. Biotechnol.* (2014), <http://dx.doi.org/10.1016/j.jbiotec.2014.08.035>
- Plotkin JB, Kudla G. **Synonymous but not the same: the causes and consequences of codon bias.** *Nat Rev Genet.* 2011;12:32–42. [Free Full Text](#)
- Shine J, Dalgarno L. **The 3'-terminal sequence of Escherichia coli 16S ribosomal RNA: complementarity to nonsense triplets and ribosome binding sites.** *Proc. Natl Acad. Sci. USA* 1974;71:1342-1346. [Free Full Text](#)
- Uebelhoer *et al.* **High-throughput, luciferase-based reverse genetics systems for identifying inhibitors of Marburg and Ebola viruses.** *Antiviral Res.* 2014 Jun;106:86-94.
- Welch M *et al.* **Design parameters to control synthetic gene expression in Escherichia coli.** *PLoS One.* 2009 Sep 14;4(9):e7002. doi: 10.1371/journal.pone.0007002.

Thank you!



- ◆ Please complete the survey
- ◆ Questions/feedback: rachel.speer@genscript.com
- ◆ Webinar Archives: www.genscript.com/webinars.html
- ◆ Request codon optimization: gene@genscript.com