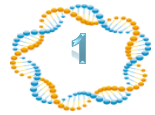


# Building DNA Libraries to Explore the Combinatorial Design Space

Dr. Yifan Li  
Senior Scientist  
GenScript USA Inc.

# Outline

---



Combinatorial Optimization for Metabolic Pathway Engineering

---



Case Studies for Combinatorial Optimization

---

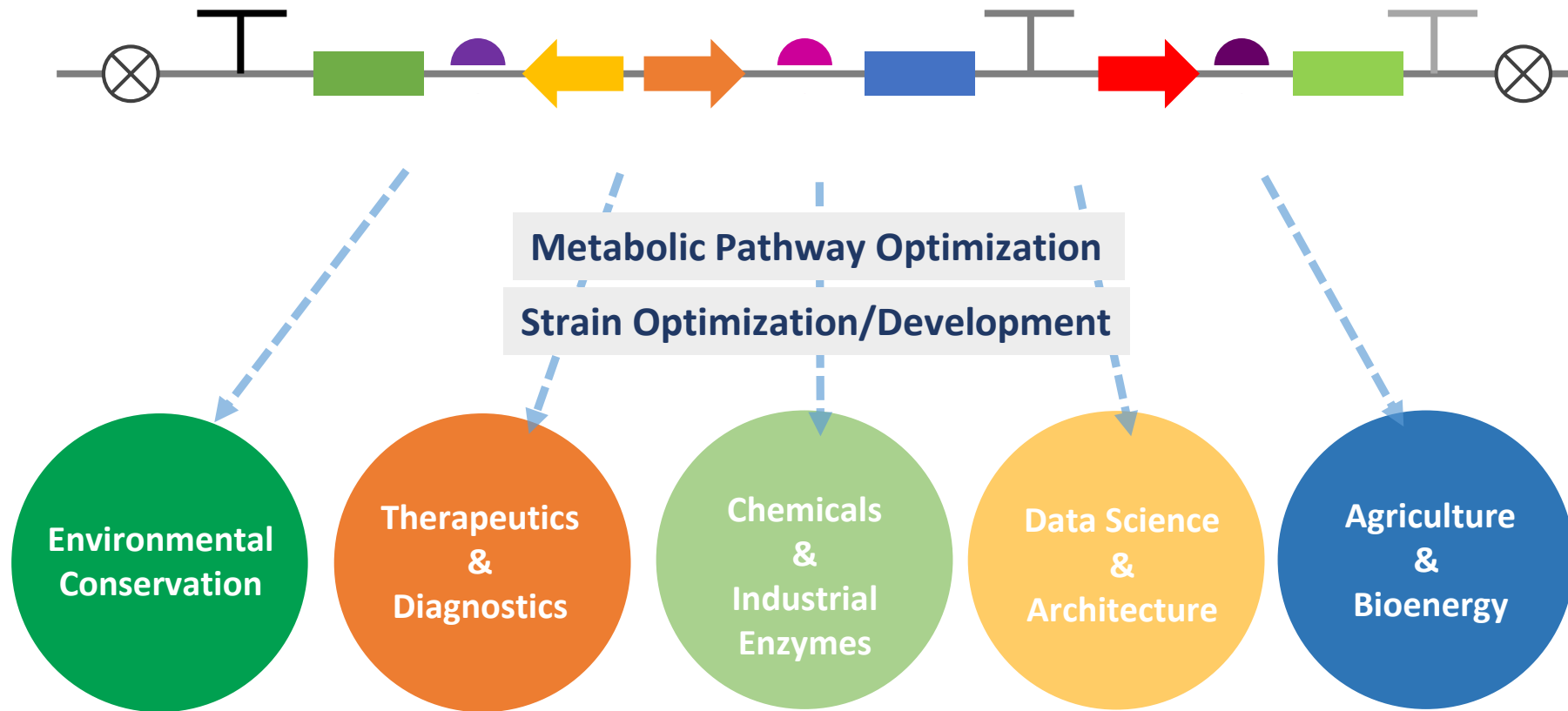


Combinatorial DNA Library Services

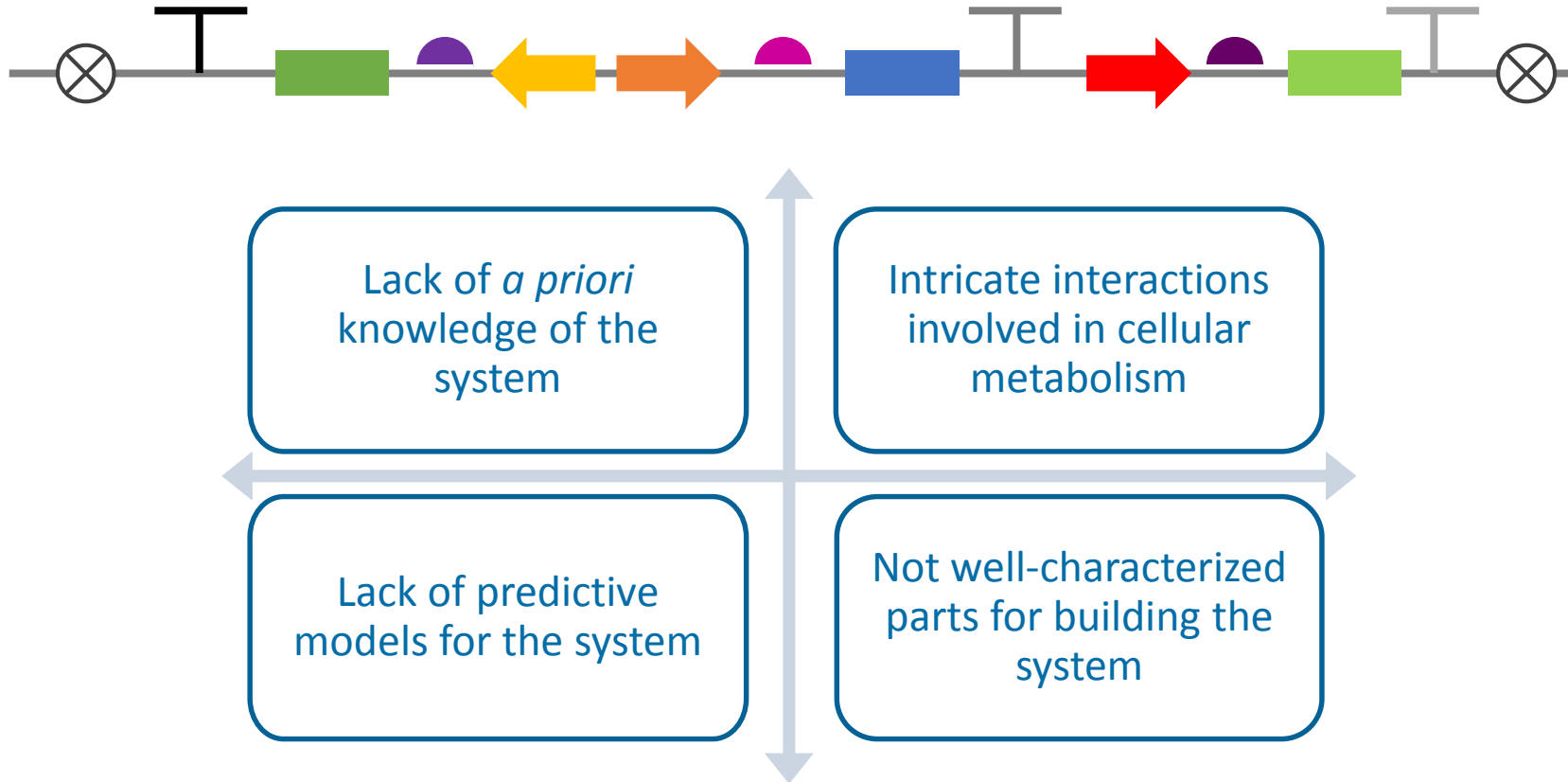
---

# Building Metabolic Pathways and Gene Circuits for Diverse Applications

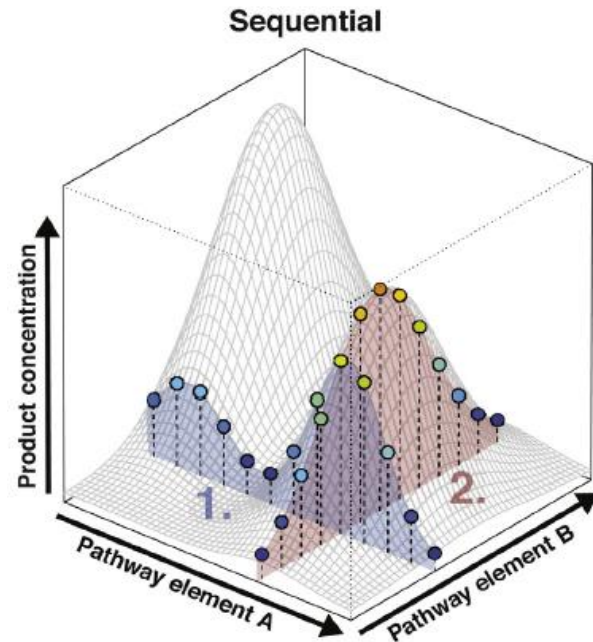
Using biological data and engineering principles to design and build novel biological systems or optimize natural systems in a more predictable and reliable manner for useful purposes.



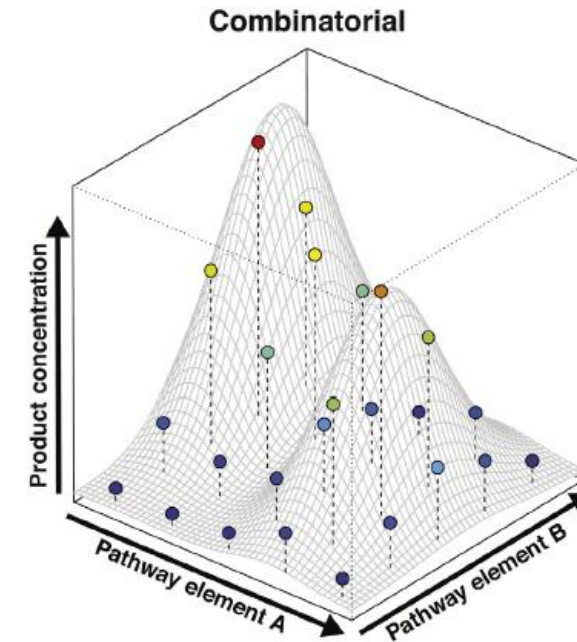
# Challenges in Engineering Biological Systems



# Sequential vs Combinatorial Optimization



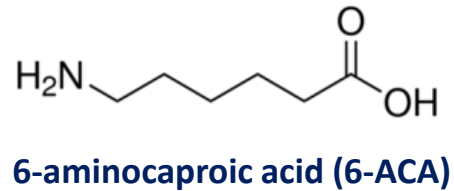
- For sequential pathway optimization, a pathway element is varied individually (blue curve) until a local optimum is identified.
- The best point is fixed as a starting point for the optimization of a second element (red curve) and so forth.



- For combinatorial pathway optimization, multiple elements are varied simultaneously.
- Combinatorial optimization allows for systematic screening of the multidimensional space and is more likely to reach a global optimum (red point).

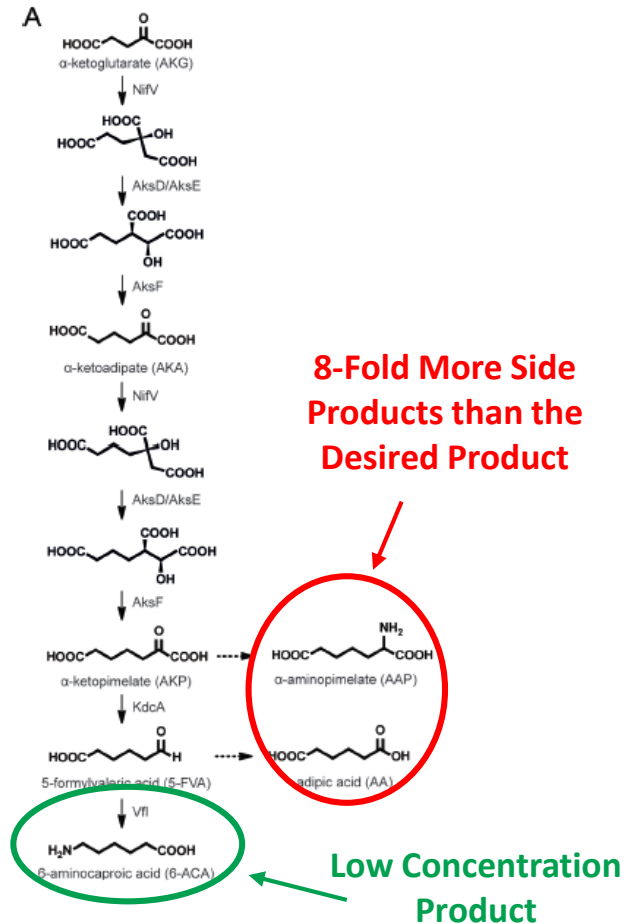
Curr Opin Biotechnol. 2017 Oct;47:142-151.

# Biosynthetic Pathway for the Production of 6-ACA



Nylon

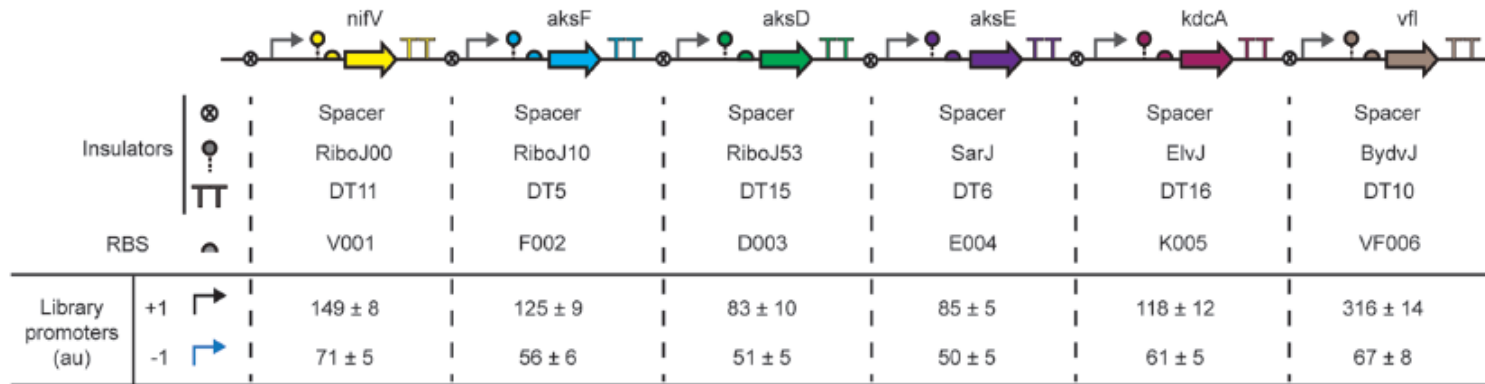
## 6-ACA Biosynthetic Pathway



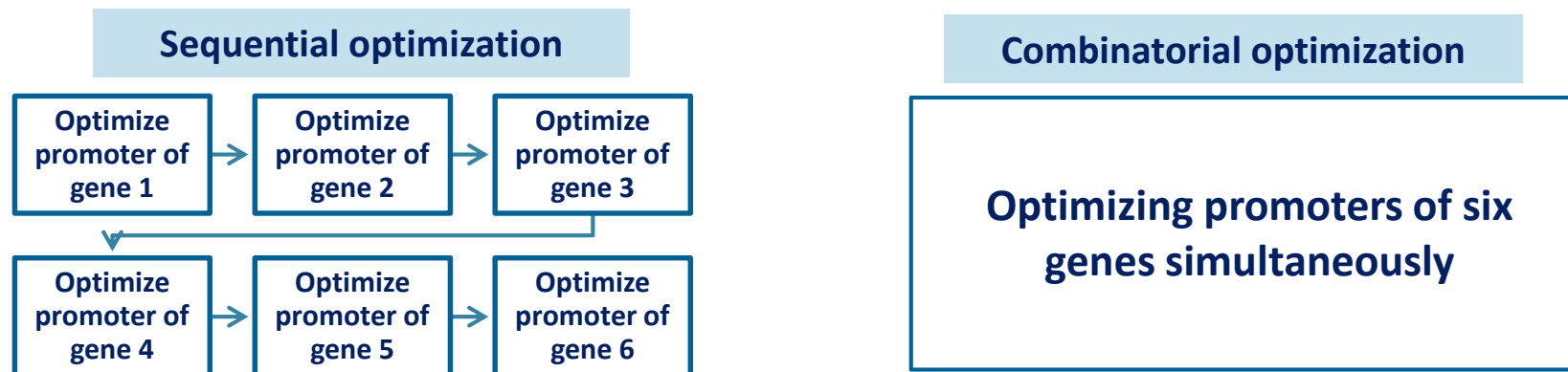
- 6-ACA is the linear form of caprolactam, which is the chemical building block of nylon-6.
- The building block is currently produced from fossil-based chemical processes, which leads to significant greenhouse gas emission.
- The natural 6-ACA biosynthetic pathway consists of six heterologous enzymes.
- The existing cell factory:
  - produced the desired product at only [8 mg/L],
  - there was 8-fold more side products than the desired product in the broth.

Nucleic Acids Res. 2015 Dec 2;43(21):10560-70.

# Pathway Optimization for 6-ACA Production



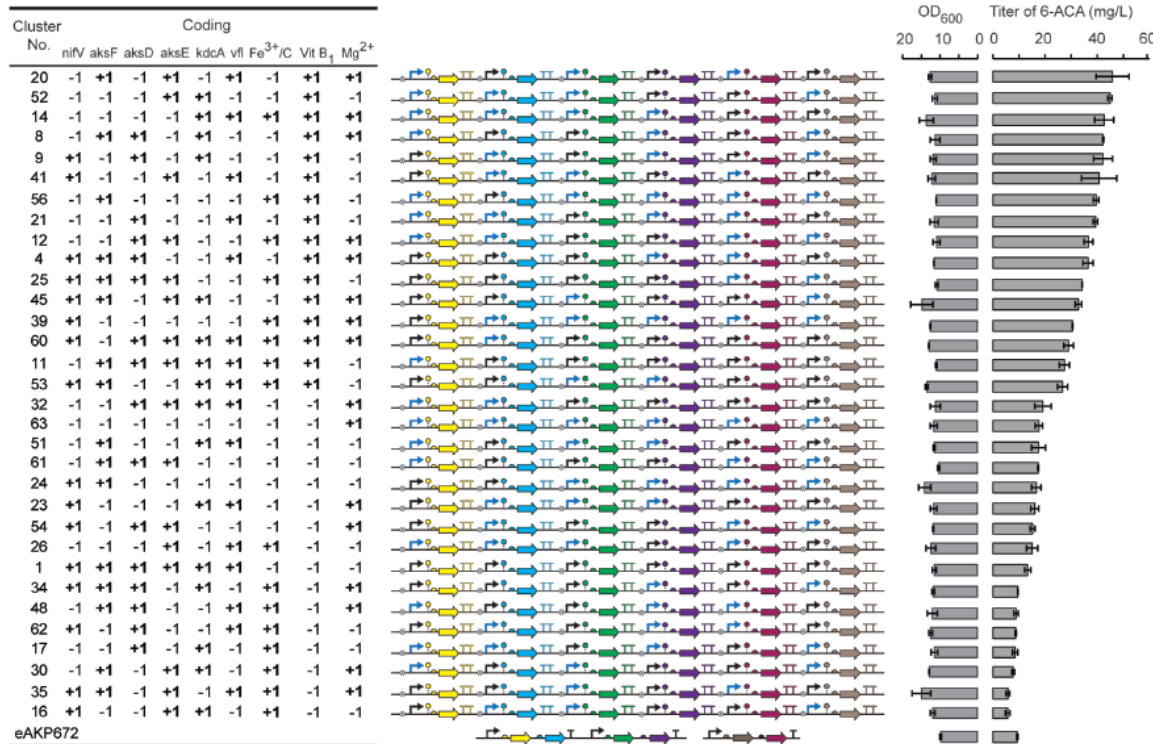
Testing strong vs weak expression for each of the six genes in the pathway



Nucleic Acids Res. 2015 Dec 2;43(21):10560-70.



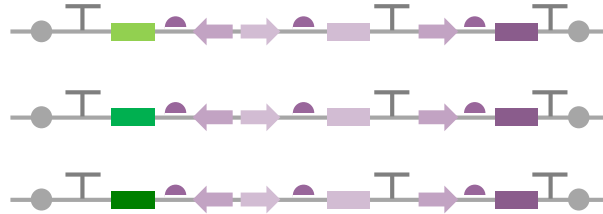
# Combinatorial Optimization for 6-ACA Production



- Three additional factors to consider in media composition:
  - [ferric citrate]
  - [vitamin B1]
  - [magnesium ion (Mg<sup>2+</sup>)]
- In total, nine factors were optimized simultaneously.
- Total library size was 2<sup>9</sup>, which equals to 512 different combinations.
- “Design of Experiment” was used to reduce combinatorial library size.
- Finally, with a single round of combinatorial optimization, 6-ACA production was increased by 5 folds.

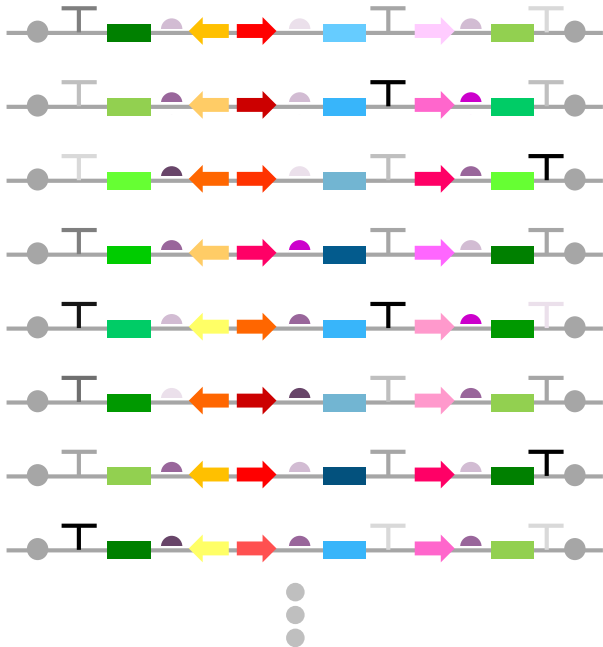


# Combinatorial Optimization Requires a Powerful Assembly Method



## Sequential optimization

- Tests:
  - less than 10 constructs at a time
  - one part at a time
- Time-consuming and costly

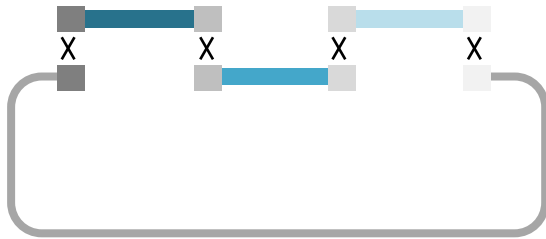


## Combinatorial optimization

- Tests:
  - hundreds and thousands constructs in parallel
  - multiple parts simultaneously
- Efficient and cost-effective

# Pathway Assembly Platforms

## *In vitro* Homologous Recombination



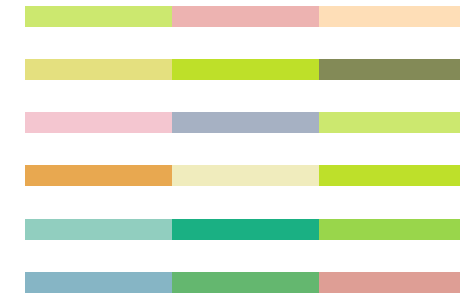
- Low throughput
- Low efficiency with >5 parts
- Expensive

## Golden Gate Assembly



- Based on Type IIs enzyme
- Has sequence limitation

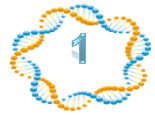
## GenScript's Platform



- High throughput
- No sequence limitation
- Seamless (no scars)
- Highly efficient with up to 10 parts and 20 kb insert size

# Outline

---



Combinatorial Optimization for Metabolic Pathway Engineering

---



Case Studies for Combinatorial Optimization

---

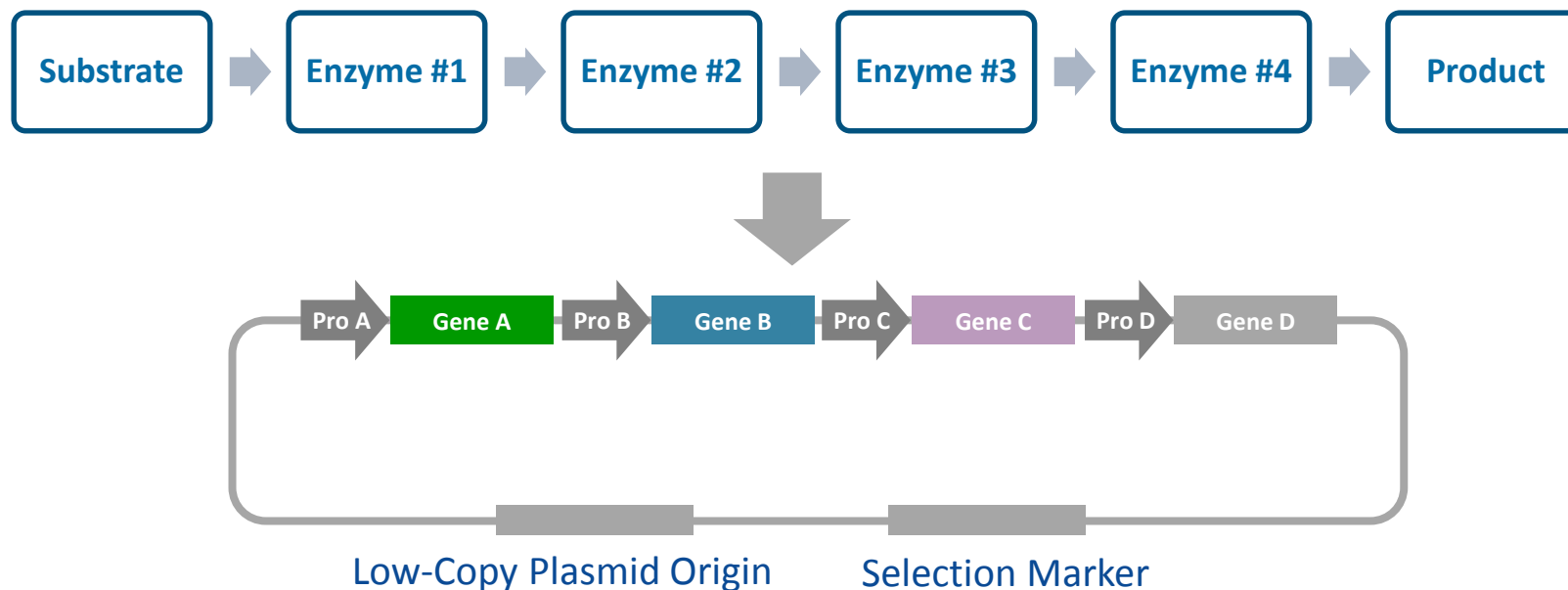


Combinatorial DNA Library Services

---

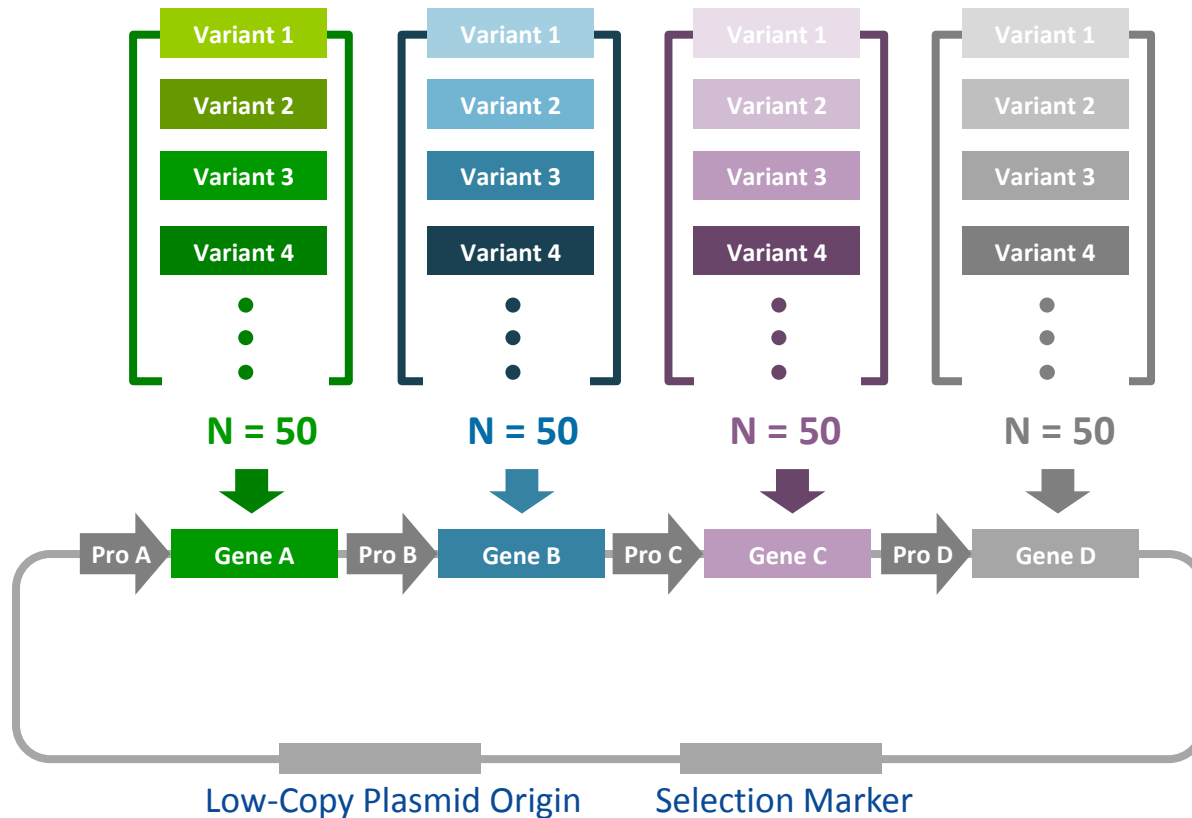
# Case #1: Metabolic Pathway Design and Construction

Objective: Improving metabolite (*i.e.* product) yield



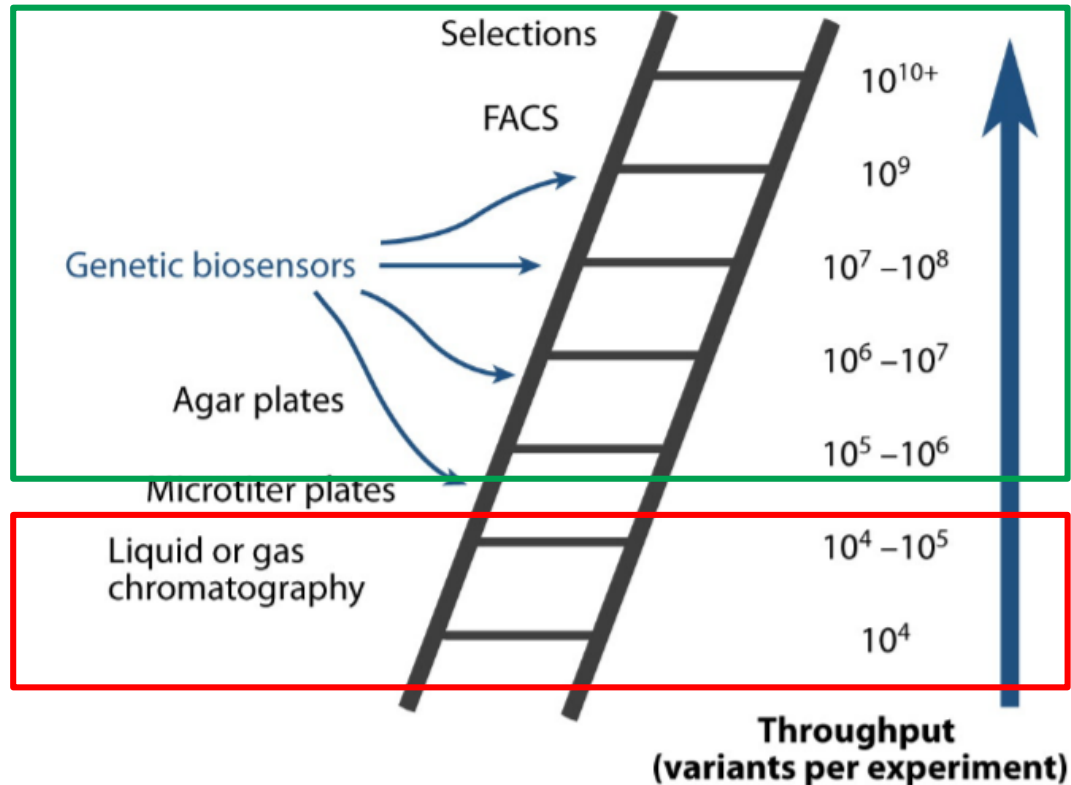
- Using yeast cells to produce the precursor for a small molecule drug
- Construct a four gene metabolic pathway in a low-copy number vector
- Using four constitutive promoters to efficiently express all four genes

# Case #1: Pathway Optimization by Testing Different Enzyme Variants



- Identified 50 variants for each pathway gene through database blasting and literature search.
- Design: Random insertion of 50 gene variants in each CDS position in the pathway
- Theoretical library size: 6,250,000

# Case #1: Throughput Screening for Different Testing Methods

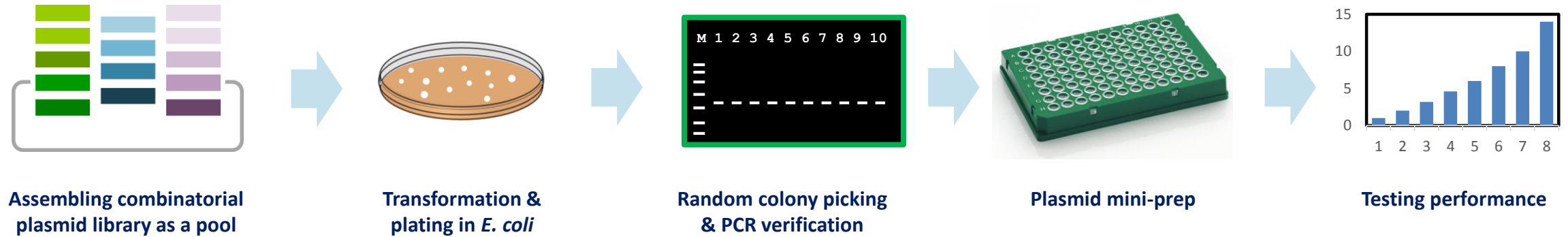


- Testing method: growth-based selection or visual selection of pigment
- Screening capacity:  $>10^6$  library size
- Explore a large library size with pooled constructs.

- Testing method: chromatography
- Screening capacity:  $<10^3$  library size
- Carefully design parameters in the project or use “Design of Experiment” to reduce library size.

Annu Rev Biochem. 2010;79:563-90.

# Case #1: Exploring Combinatorial Library by Randomly Picking Clones



## Workflow

- Library construction as a pool – putting all fragments into the reaction
- Randomly tested 2,000 colonies; ~1600 of them have inserts, indicating an 80% positive rate.
- Mini-prepped all plasmids with HT platform, followed by transformation of plasmids into appropriate expression host to test performance.

## Key Points:

- By not assembling and sequencing all combinations, lots of money and effort was saved.
- By testing a sub-library, enough time was saved for HPLC testing.
- Colony PCR verification ensures that every tested clone was positive (*i.e.* contained the designed components).



# Case #1: Testing Library Diversity to Ensure Maximum Coverage

Combination #	Enzyme A#	Enzyme B#	Enzyme C#	Enzyme D#
1	45	25	40	19
2	8	24	16	30
3	49	37	32	6
4	9	41	31	30
5	36	10	22	8
6	44	13	37	4
7	15	44	14	2
8	23	28	4	13
9	15	44	12	20
10	13	12	29	47
11	32	37	3	40
12	36	18	31	5
13	29	1	3	17
14	21	23	27	22
15	2	28	36	25
16	38	27	35	45
17	4	28	50	35
18	40	32	35	35
19	4	16	18	9
20	18	8	39	4
21	33	10	28	7
22	48	40	40	44
23	31	34	16	37
24	21	6	22	9

Combination #	EnzymeA #	EnzymeB #	EnzymeC #	Enzyme D#
25	43	1	6	26
26	14	36	48	9
27	10	25	26	15
28	34	24	10	4
29	32	1	32	50
30	11	30	21	7
31	2	43	46	28
32	38	46	1	49
33	13	18	47	3
34	37	50	40	5
35	8	17	42	41
36	5	22	49	38
37	5	19	1	10
38	20	4	48	38
39	47	27	31	49
40	14	46	23	2
41	16	47	24	49
42	13	29	14	11
43	36	37	19	23
44	3	32	2	36
45	45	34	2	37
46	35	5	3	14
47	35	22	15	19
48	15	36	39	49

Analyzing the sequence of 48 randomly-picked clones demonstrated that 100% diversity was achieved.

# Case #1: Screening for the Best Clones

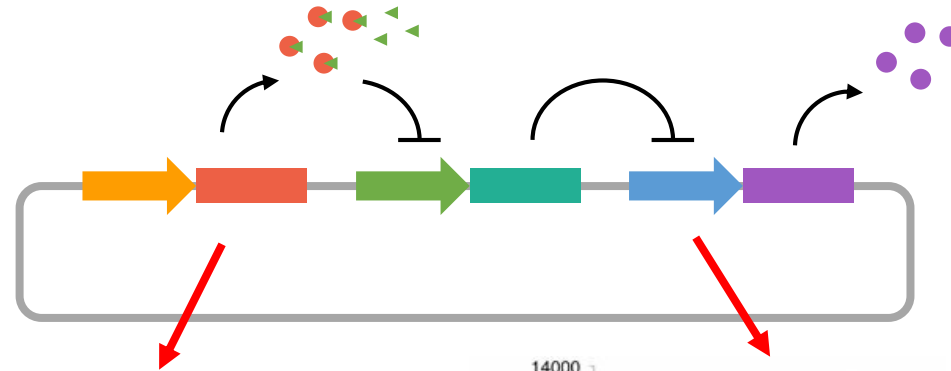
## Top 24 Variants, Ranked by Production Level

Combination #	Enzyme A#	Enzyme B#	Enzyme C#	Enzyme D#	Production (Arbitrary unit)
1	13	49	11	23	
2	13	49	5	44	
3	10	46	11	17	
4	37	46	11	7	
5	26	42	11	44	
6	13	33	5	45	
7	44	44	27	40	
8	27	16	5	31	
9	14	46	38	9	
10	38	28	5	44	
11	49	42	30	46	
12	14	49	27	14	
13	37	16	5	6	
14	16	6	12	19	
15	13	49	11	45	
16	37	42	1	6	
17	32	36	39	23	
18	26	6	27	12	
19	37	42	4	20	
20	35	49	5	33	
21	42	17	22	31	
22	15	16	2	7	
23	3	34	3	8	
24	32	28	11	12	

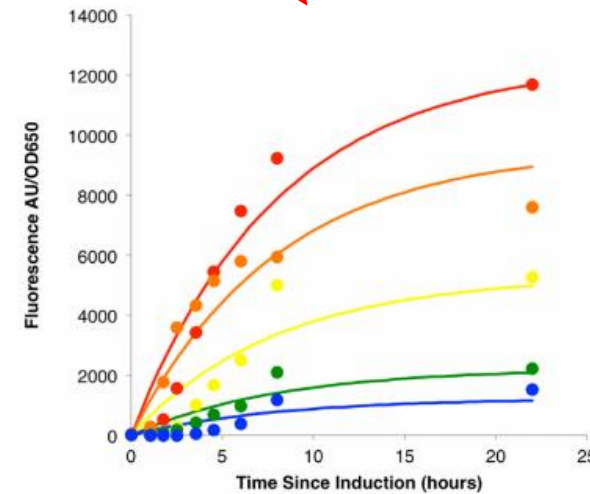
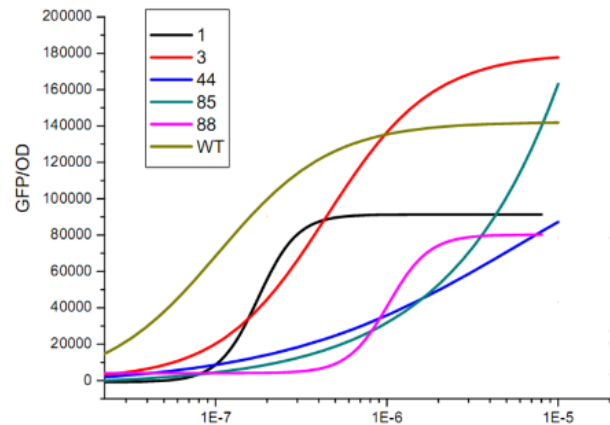
- Tested all 2000 randomly-picked clones with HPLC analysis.
- Sequence analyzed the top 96 variants. (Top 24 variants are listed here.)
- Some of the enzyme variants were highly enriched in the good-performing clones.
- These good-performing enzyme variants can be further combinatorically tested in the next round of optimization.

## Case #2: Optimizing Gene Circuits for Therapeutic Applications

Objective: Production of a functional protein upon sensing specific environmental signals

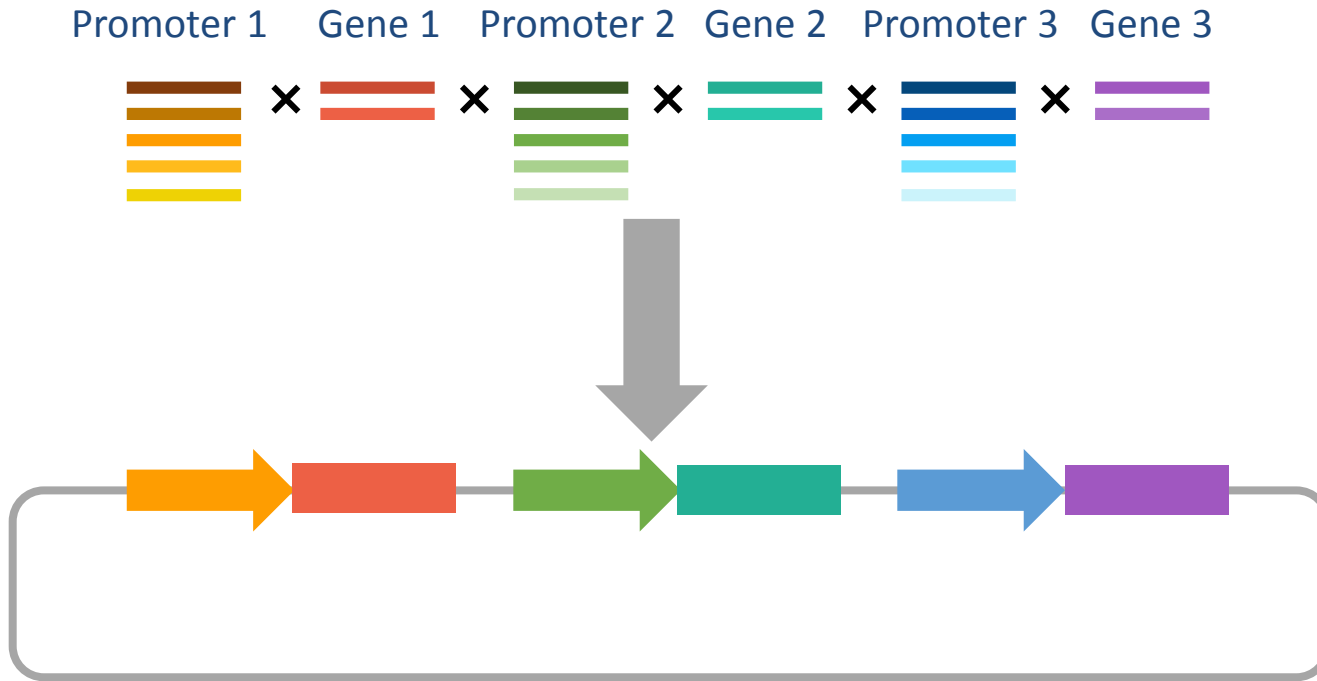


Different dose response curves for different regulators was needed.



Expression intensity could affect performance of the circuit.

## Case #2: Optimization of Gene Circuits



### Project scope

- Testing five promoters with strength levels from very weak to very strong for gene expression
- Testing different versions of sensors and regulators
- Theoretical combination number:  
 $5 \times 2 \times 5 \times 2 \times 5 \times 2 = 1,000$

# Case #2: Construction & Testing of Each Combination Individually



## Workflow:

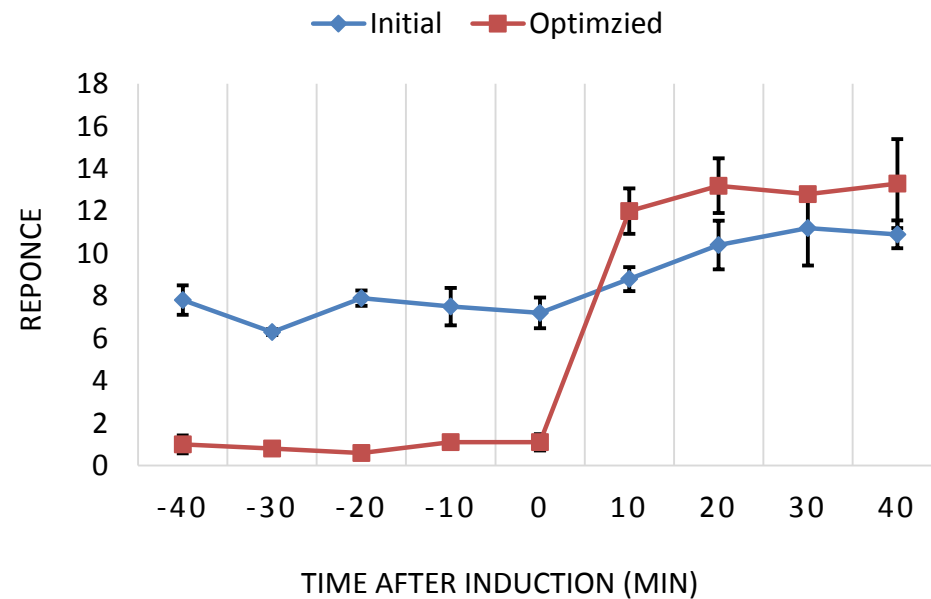
- Assembled 1,000 construct individually and acquired 943 perfect clone in the first run.
- Constructed the remaining 57 clones in the second run.

## Key Points:

- Individual building and testing of every construct allowed for direct evaluation of each construct.
- Suitable for thoroughly explore the combinatorial design space
- Compatible with “Design of Experiment”

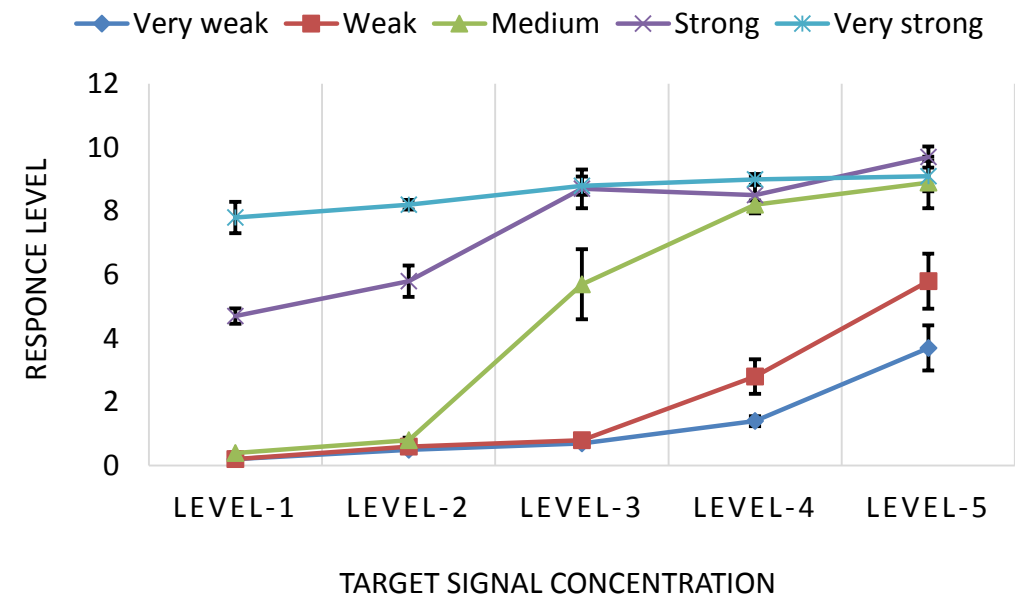
## Case #2: Optimized Circuit exhibited Improved Performance

Response Curve after Induction



The optimized circuit exhibited stronger and more prompt response behavior than the initial.

Inducer Dose Response for Promoter-3 Strength



The medium-strength promoter-3 showed the best dose-response behavior.

# Summary of Case Studies

## Case #1

### Representative Combinatorial DNA Library

---

- Library construction as a pool
- Random colony-picking for downstream testing
- Saves a lot of money and effort
- Suited for the initial phase of biological optimization

## Case #2

### Arrayed Combinatorial DNA Library

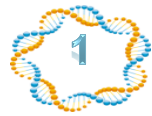
---

- Building and testing of every construct individually
- Suitable for thoroughly exploring the combinatorial design space
- Specifically suitable for optimizing non-linear systems
- Compatible with “Design of Experiment”



# Outline

---



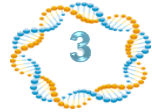
Combinatorial Optimization for Metabolic Pathway Engineering

---



Case Studies for Combinatorial Optimization

---



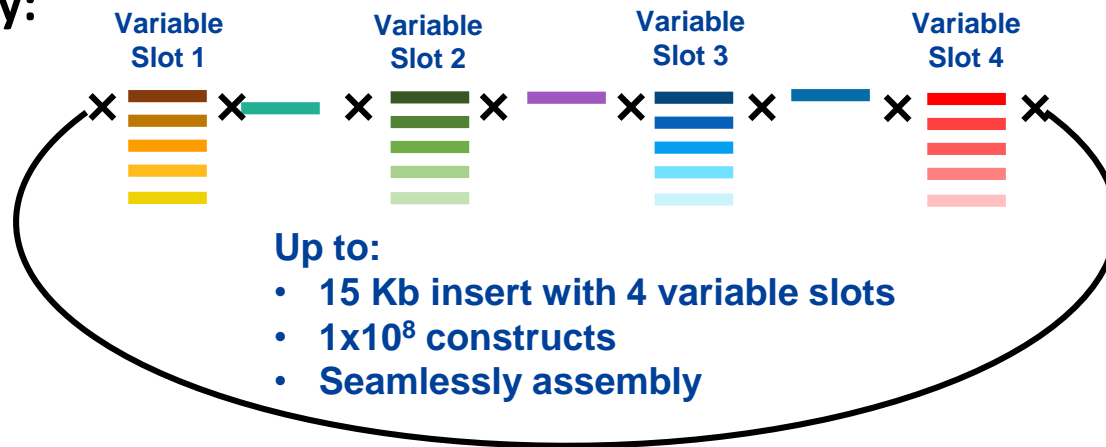
Combinatorial DNA Library Services

---

# Combinatorial DNA Libraries

Providing a **highly customizable** approach for **accelerating** the **build phase** of your **metabolic pathway** and **microbial strain engineering process**!

## Our Capability:



## Applications:

- *Metabolic Pathway & Microbial Strain Engineering*
- *Gene Expression Regulation*
- *Protein & Antibody Engineering*

# Combinatorial DNA Library Services

## Pooled Combinatorial DNA Library



### Ideal for

- $> 10^4$  throughput screening
- For screening platforms that are not sensitive to the presence of negative clones

### Service Features

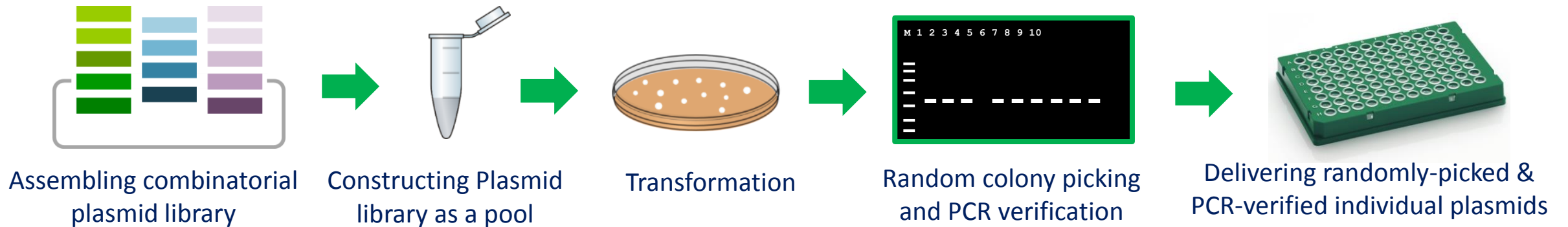
- Cloned into pUC57 or custom vector with 4  $\mu$ g in quantity
- Delivering pooled a plasmid library with up to  $1 \times 10^8$  library size

### QC Standards

- PCR verification of more than 48 clones to determine positive rate
- Sequence verification of 24 positive clones with a guarantee on more than 85% diversity

# Combinatorial DNA Libraries

## Representative Combinatorial DNA Library



### Ideal for

- $10^2$  -  $10^4$  throughput screening
- Screening a pool with no concern for the presence of exact sequence in each clone

### Service Features

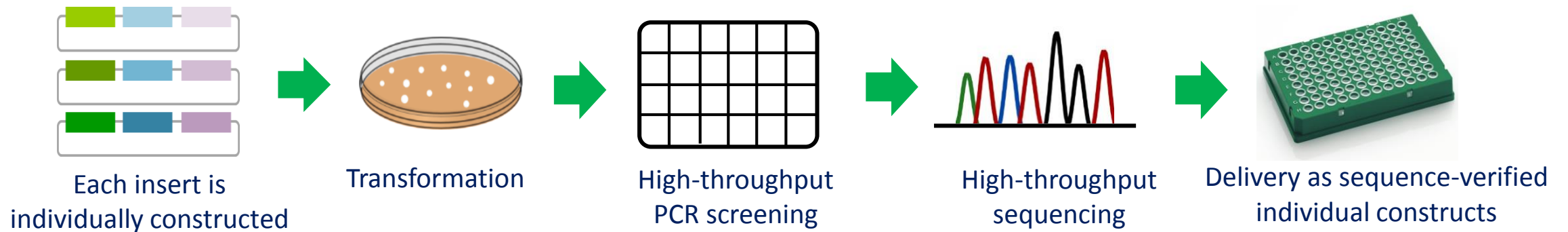
- Guaranteeing that every delivered construct contains all designed parts or modules
- Delivering up to 10,000 randomly-picked and PCR-verified individual constructs
- Cloned into pUC57 or custom vector with 4  $\mu$ g in quantity

### QC Standards

- PCR verification of all delivered plasmids
- Sequence verification of 24 positive clones with a guarantee on more than 85% diversity

# Combinatorial DNA Libraries

## Arrayed Combinatorial DNA Library



**Ideal for** Testing every single design in your library

### Service Features

- Enabling the design of every construct in the library in any combination
- Delivering up to 10,000 sequence-verified individual constructs ready for transformation
- Cloned into pUC57 or custom vector with 4 µg in quantity

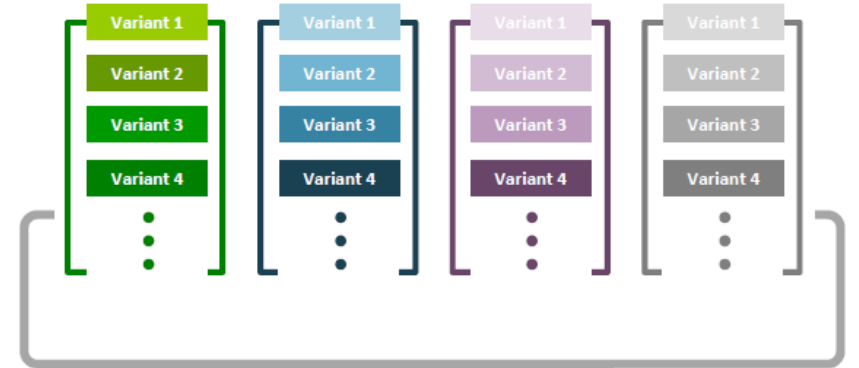
### QC Standards

Each delivered plasmid is sequence-verified.

# Combinatorial DNA Libraries: Advantages

*Let DNA Building Experts Speed Up Your Metabolic Pathway  
and Microbial Strain Engineering Process!*

- ✓ **One-stop, high-throughput** solution
- ✓ **Highly-customizable**
  - In pool or individual formats
  - Up to  $1 \times 10^8$  constructs (pooled library)
  - Up to 4 variable slots for 15 kb inserts
- ✓ **Faster and more economical** compared to your in-house operations
- ✓ **Seamlessly assembled** with advanced methods
- ✓ **Expert advising** on all project plans by our Ph.D.-level scientists



# Thank you!

For any questions,  
please visit:

[www.genscript.com/combinatorial-DNA-library.html](http://www.genscript.com/combinatorial-DNA-library.html)

or email:

[clairez@genscript.com](mailto:clairez@genscript.com)  
[gene@genscript.com](mailto:gene@genscript.com)



Gene



Peptide



Protein



Antibody



Discovery



Catalog Products