

# A Guide to DNA Assembly for Drug Discovery

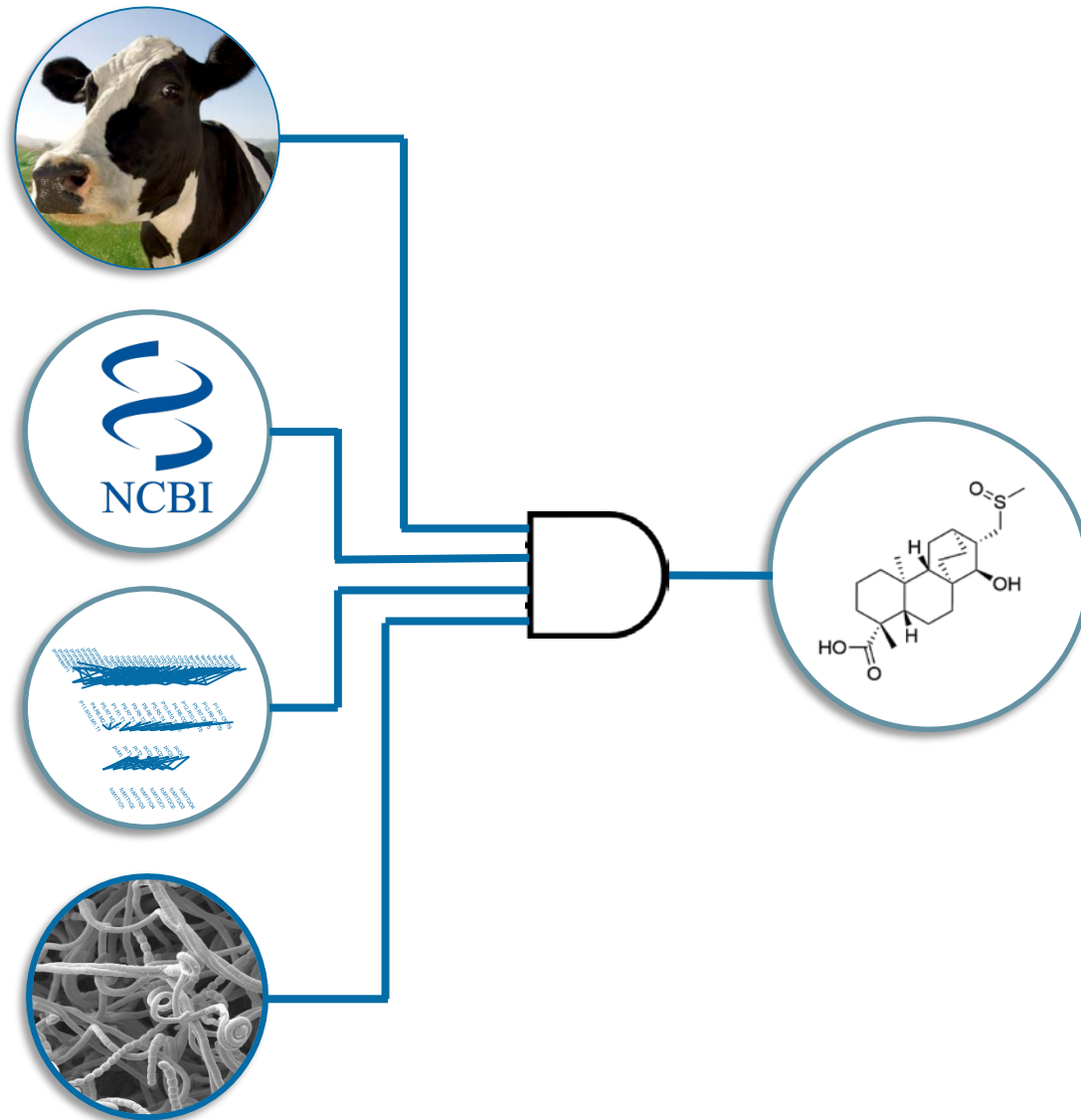
Michael Smanski, Ph.D.  
Assistant Professor  
University of Minnesota

May 15, 2018

# Leveraging DNA Assembly for Natural Product Biosynthesis

Mike Smanski  
University of Minnesota

May 15, 2018



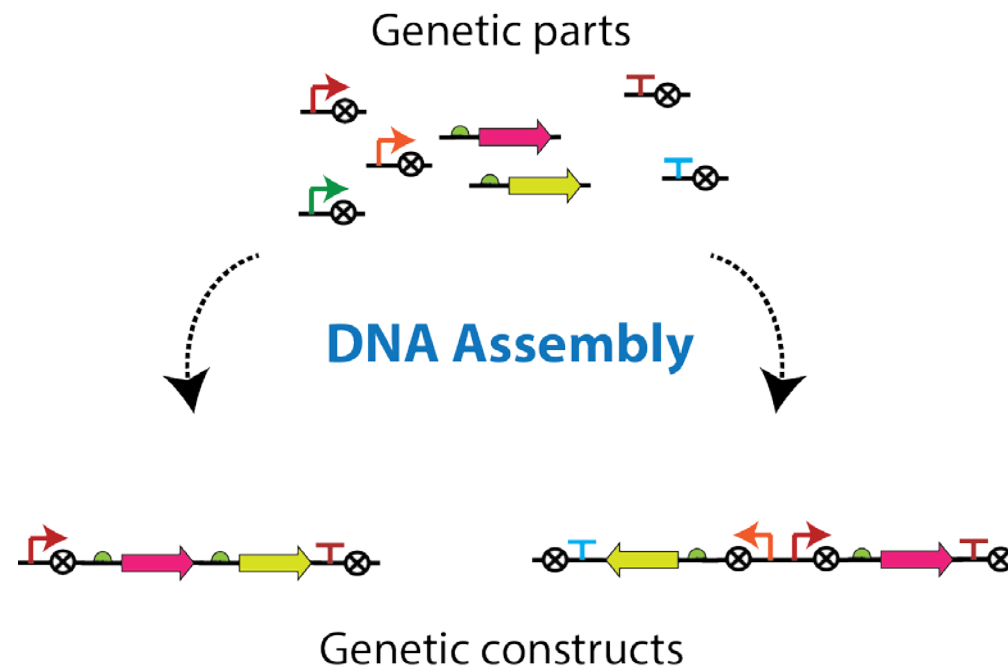
# Goals for the Webinar

- Introduction to **key concepts** in DNA assembly as it relates to engineering metabolic pathways
- Argument for the **importance of combinatorial DNA assembly** projects in natural product biosynthesis
- **Example application** to illustrate how DNA assembly methods were used to produce a valuable natural product scaffold
- **New directions/questions** that are important for pushing genetic engineering into more complex multi-part systems

# “DNA Synthesis” versus “DNA Assembly”

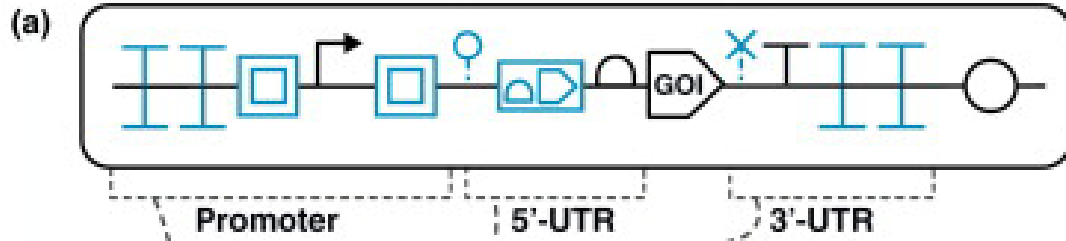


**DNA Synthesis:** de novo construction of oligonucleotides and larger molecules from nucleotide monomers



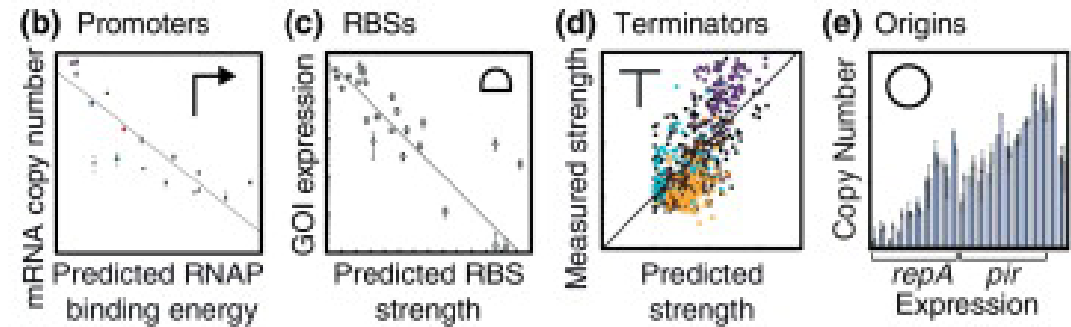
**DNA Assembly:** combinatorial concatenation of pre-synthesized parts to produce functional genetic constructs

# What do We Mean by “Genetic Parts”?

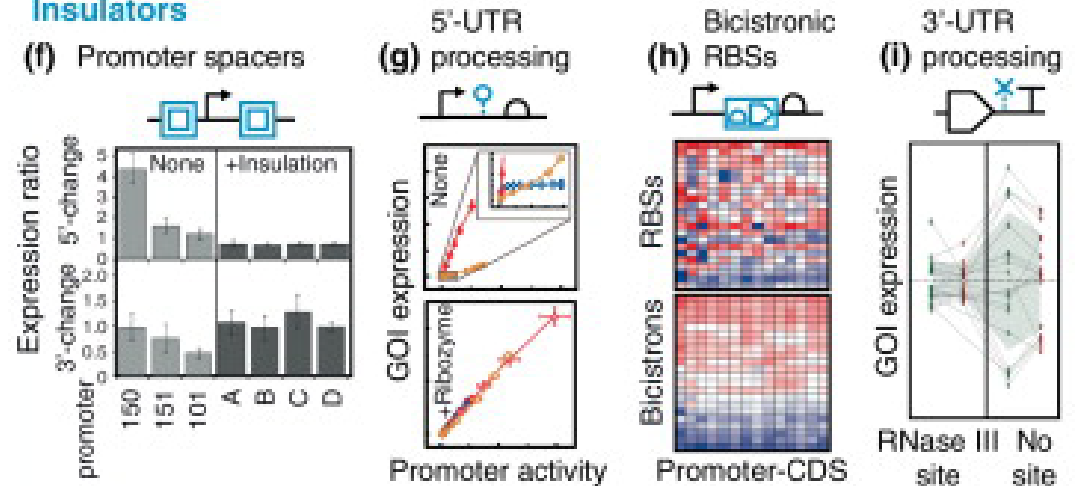


**Genetic parts:** a sequence of DNA that encodes a biological function or behavior. This can include information storage (e.g. CDS), transcription or translation control (e.g. promoter), etc. Genetic parts can be combined to make more complex functional units.

## Tuning Knobs



## Insulators

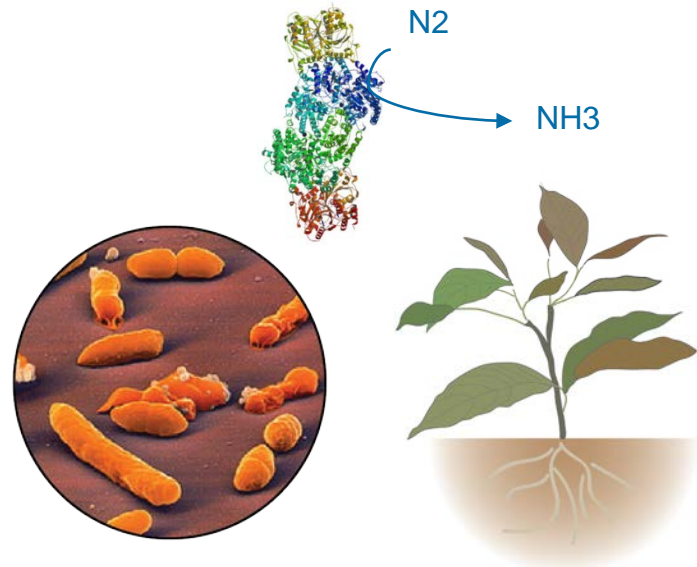


Nielsen AAK, Segall-Shapiro TH, Voigt CA *Curr. Opin. Chem. Biol.* (2013)

# The iGEM Registry of Standard Biological Parts

The screenshot shows the iGEM Registry website interface. At the top, there is a navigation bar with links for 'iGEM', 'wiki tools', 'search', 'PRODUCTION 2017 SERVER', and a 'login' button. The main heading is 'Registry of Standard Biological Parts'. Below this is a secondary navigation bar with links for 'tools', 'catalog', 'repository', 'assembly', 'protocols', 'help', and 'search', along with a search input field containing 'BBa\_'. The central content area features a large image of stacked cardboard boxes with the iGEM logo. A red banner is overlaid on the image with the text '2018 Distribution is shipping!'. To the right of the image is a 'Registry News' sidebar with a list of links: 'Registry Release', 'Registry 6.0', 'Report Bugs', 'Request Features', 'News Archive', and 'Feature Box Archive'. Below the sidebar is an 'Other' section with links for 'Registry API', 'Safety', and 'Videos'. At the bottom of the page, there are three boxes: 'iGEM 2018 Team Registration' (with subtext 'Late registration for iGEM 2018 is now...'), '2018 Distribution Kits' (with subtext 'The 2018 Distribution Kits have started...'), and 'Registry Updates' (with subtext 'The Registry will be undergoing updates...').

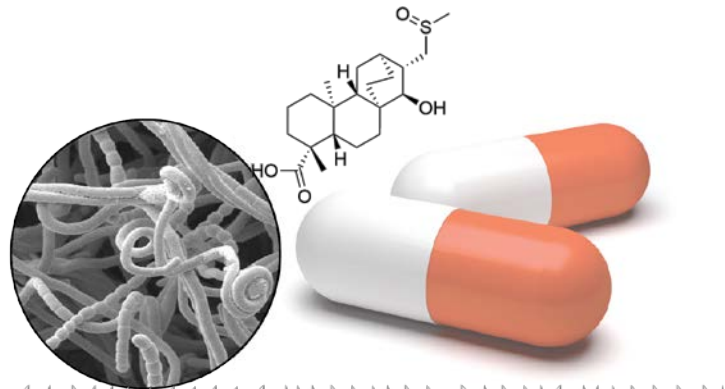
# Important Biological Capabilities Require Massively Multi-Part Genetic Systems



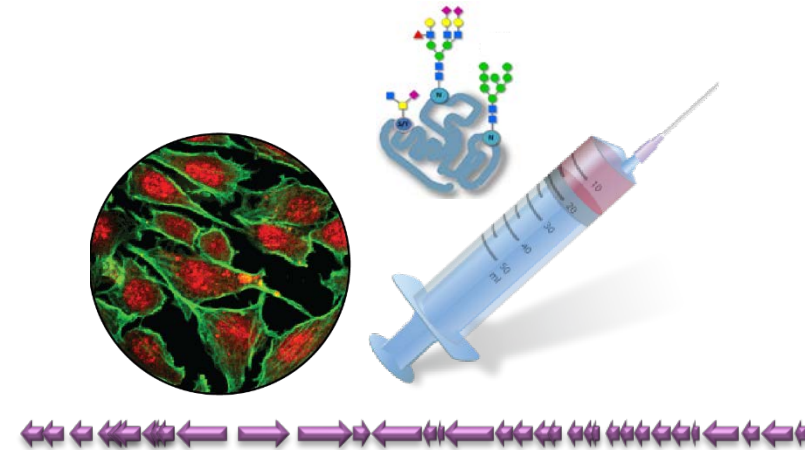
**Nitrogen Fixation**



**Electron transfer**

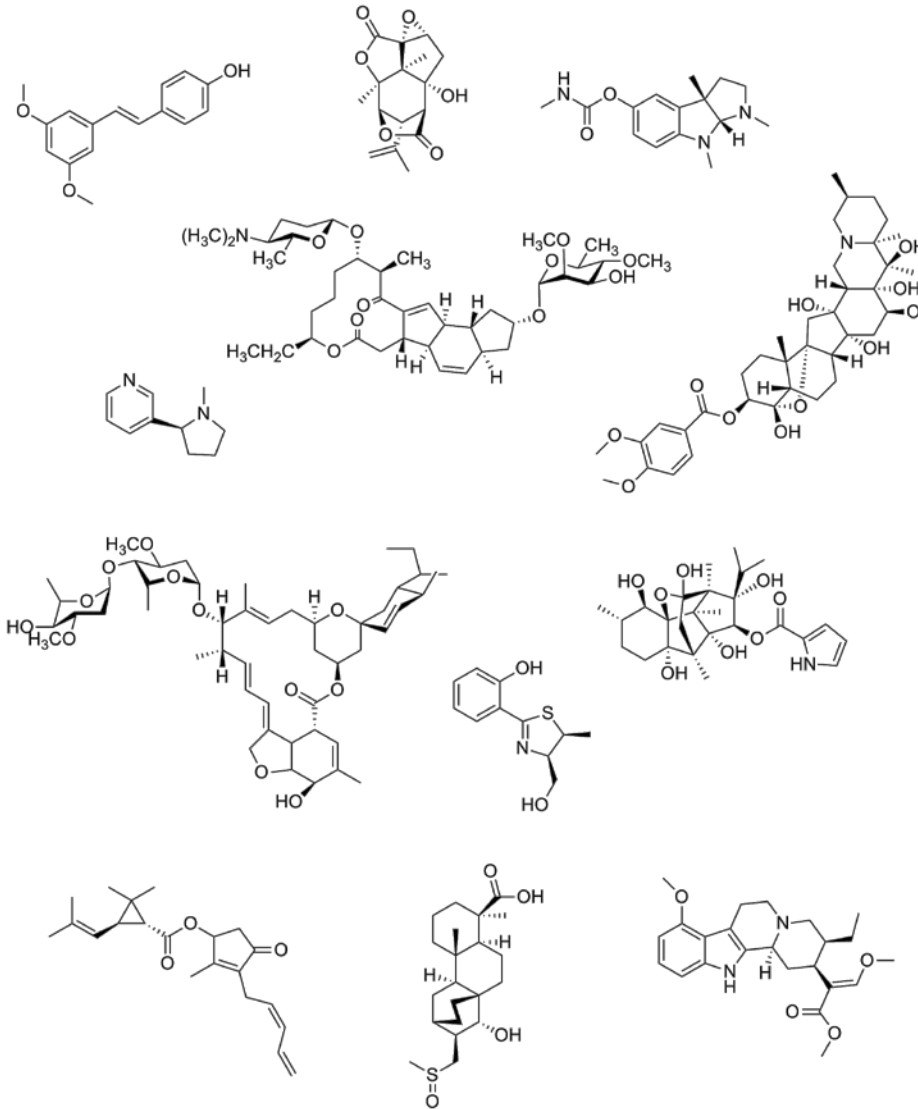


**Natural product biosynthesis**



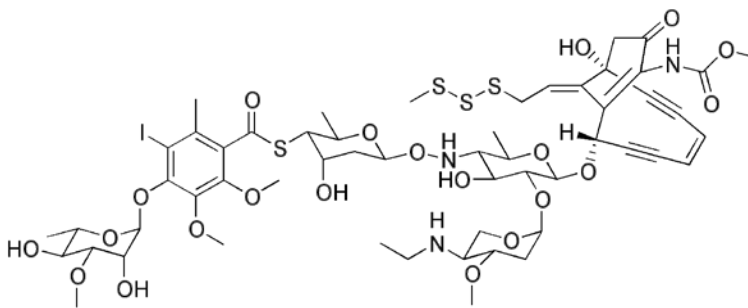
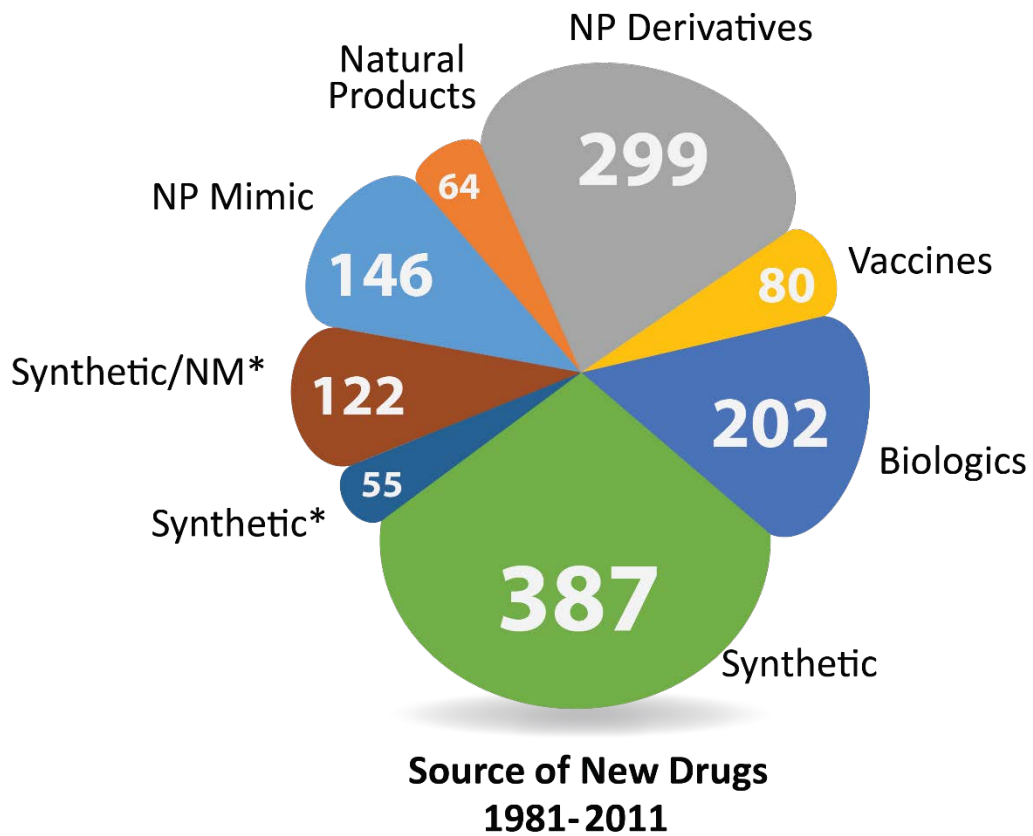
**Protein glycosylation**

# What are “Natural Products”

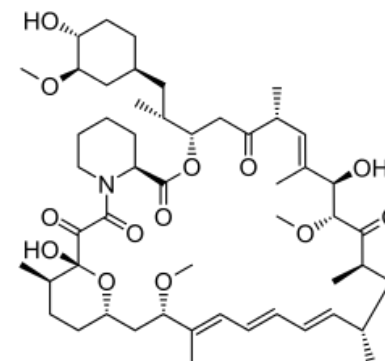




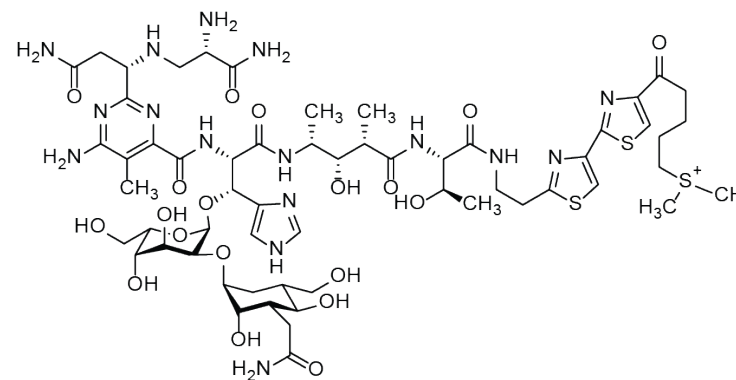
# Natural Products are Privileged Structures for Drug Discovery



Calicheamicin (Mylotarg®)



Rapamycin (Rapamune®)



Bleomycin (Blenoxane®)

Newman & Cragg (2012) *J. Nat. Prod.* **75**:311-335

# Heterologous Production of Natural Products

Native host

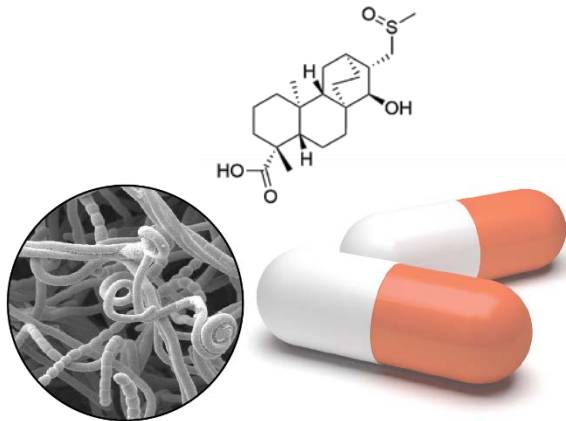
*S. platensis*



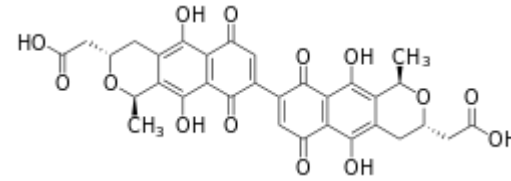
Direct gene cluster transfer

Heterologous host

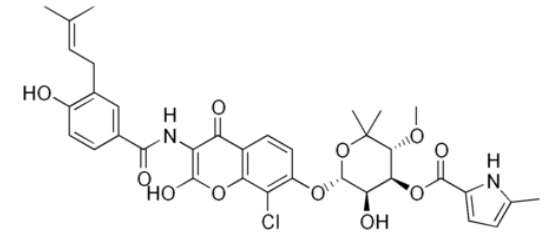
*S. lividans*



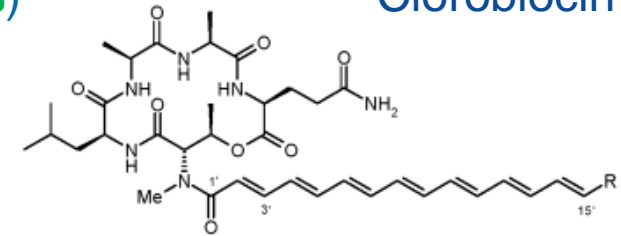
Natural product biosynthesis



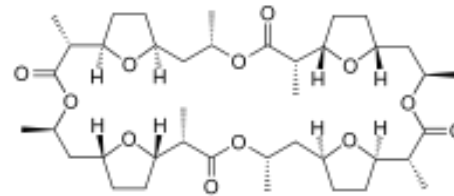
Actinorhodin (100%)



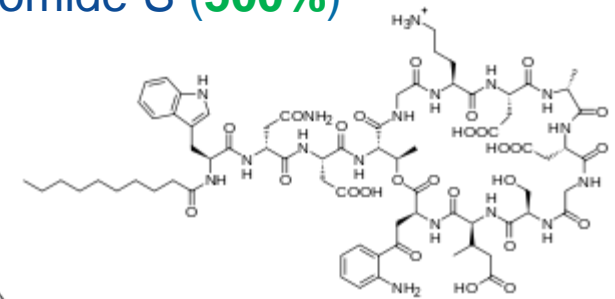
Clorobiocin (104%)



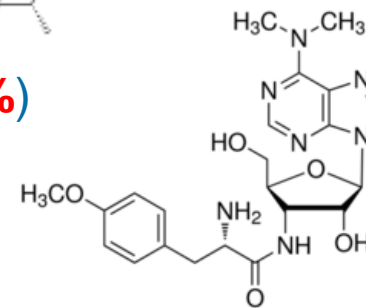
Myxochromide S (500%)



Nonactin (25%)



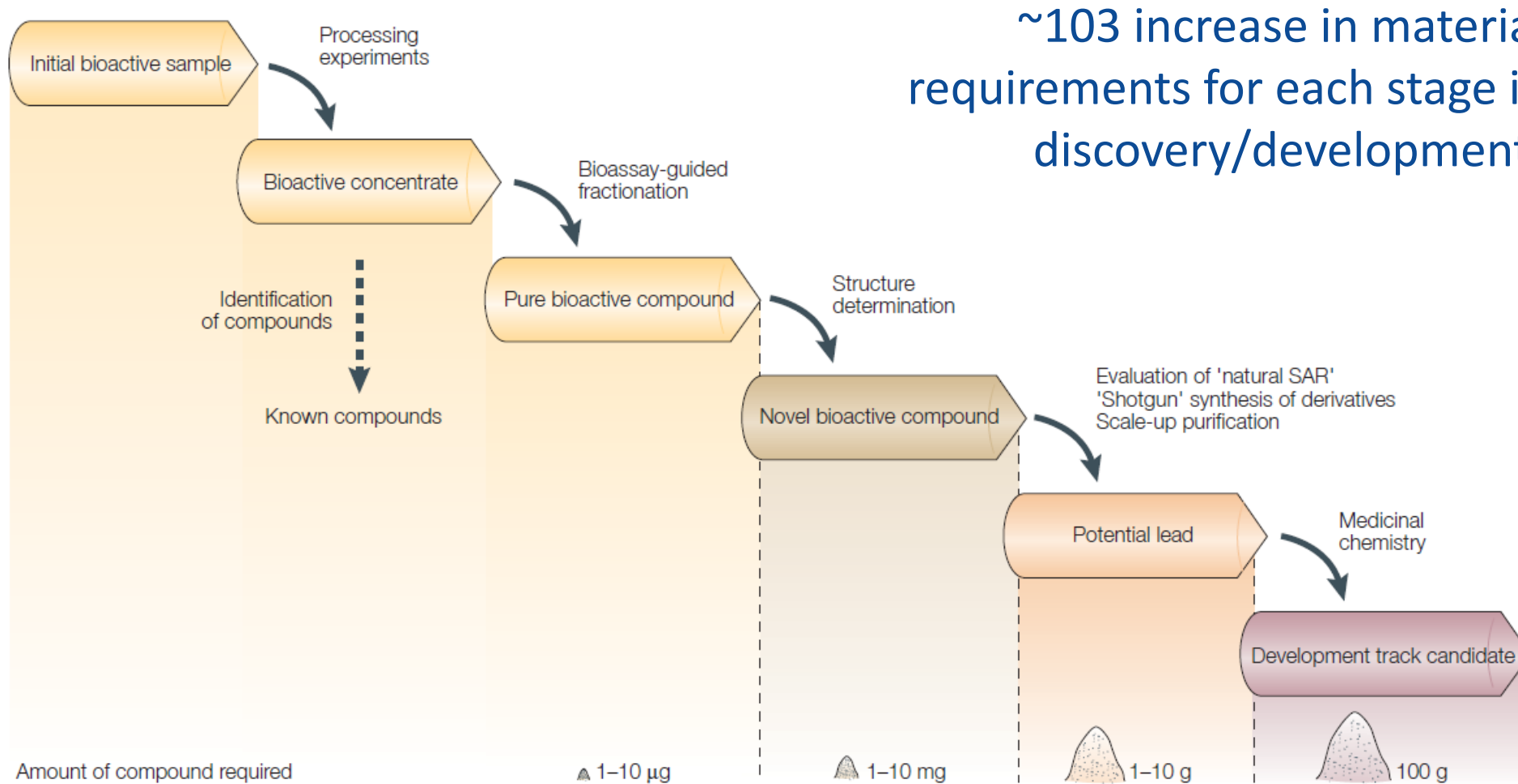
Daptomycin (2%)



Puromycin (10%)

Expert Opin. Drug Discov. (2006) 1(5)

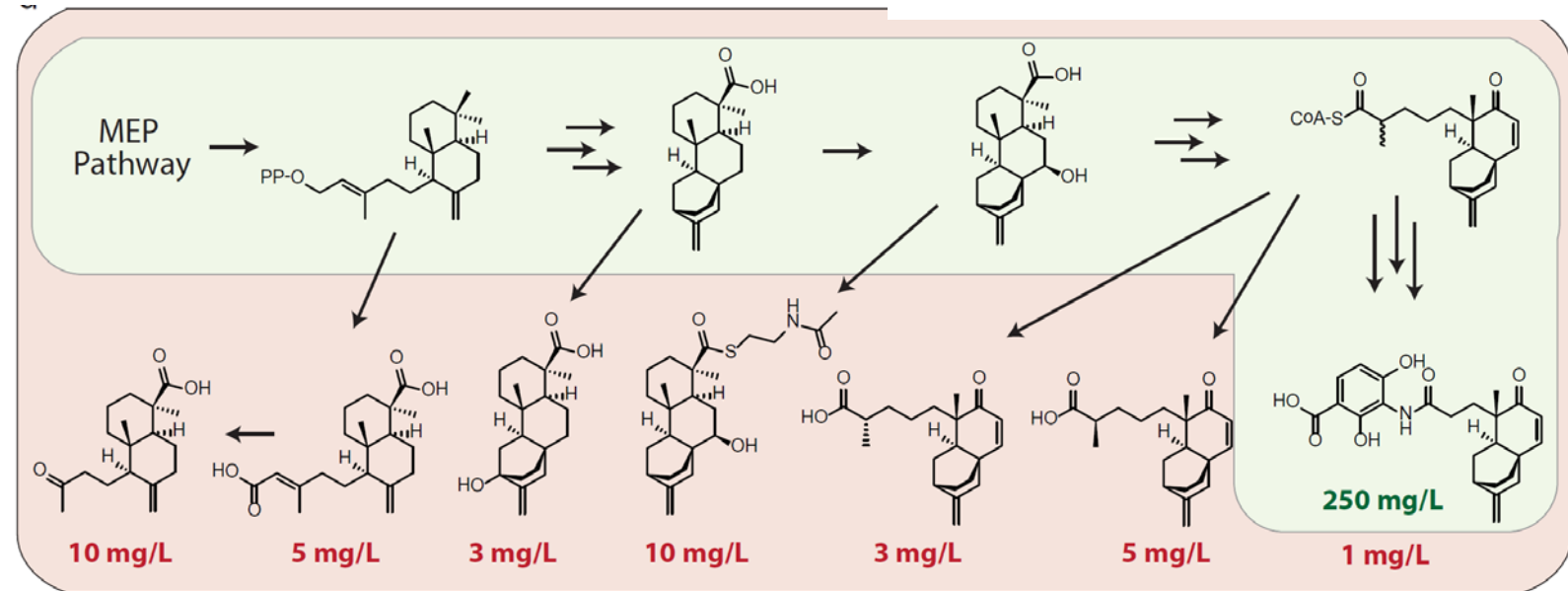
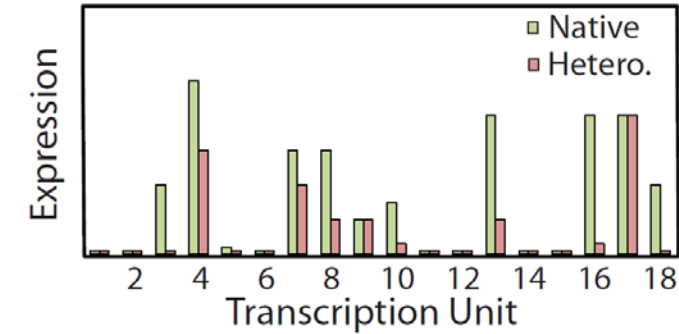
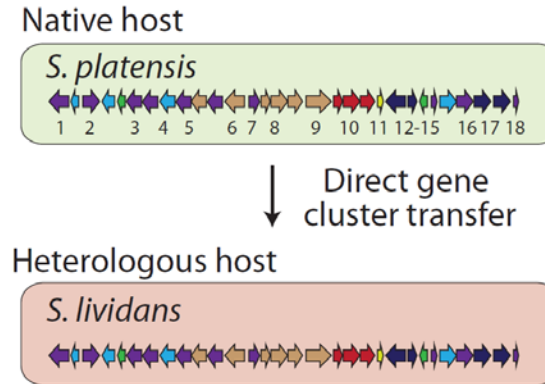
# Drug Discovery Pipelines Require Access to Hundreds of Grams of Material



Koehn and Carter (2005) *Nat. Rev. Drug Disc.* 4:206-220

# Gene Expression is Finicky but Important for Heterologous Production

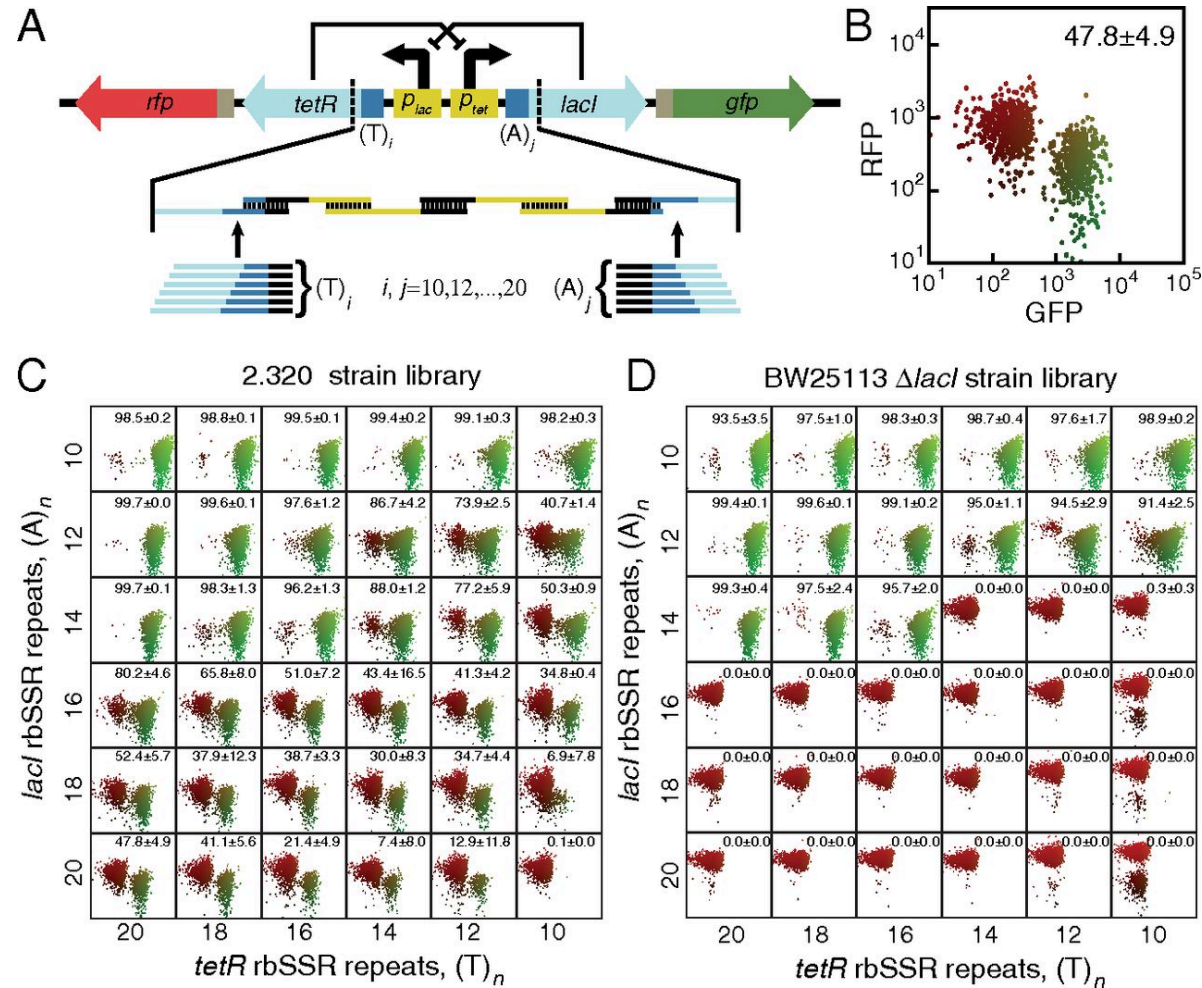
- Gene expression changes upon host transfer



Smanski et al. Shen (2012) *J. Nat. Prod.* **75**:2158-2167

# Gene Expression is Finicky but Important for Heterologous Production

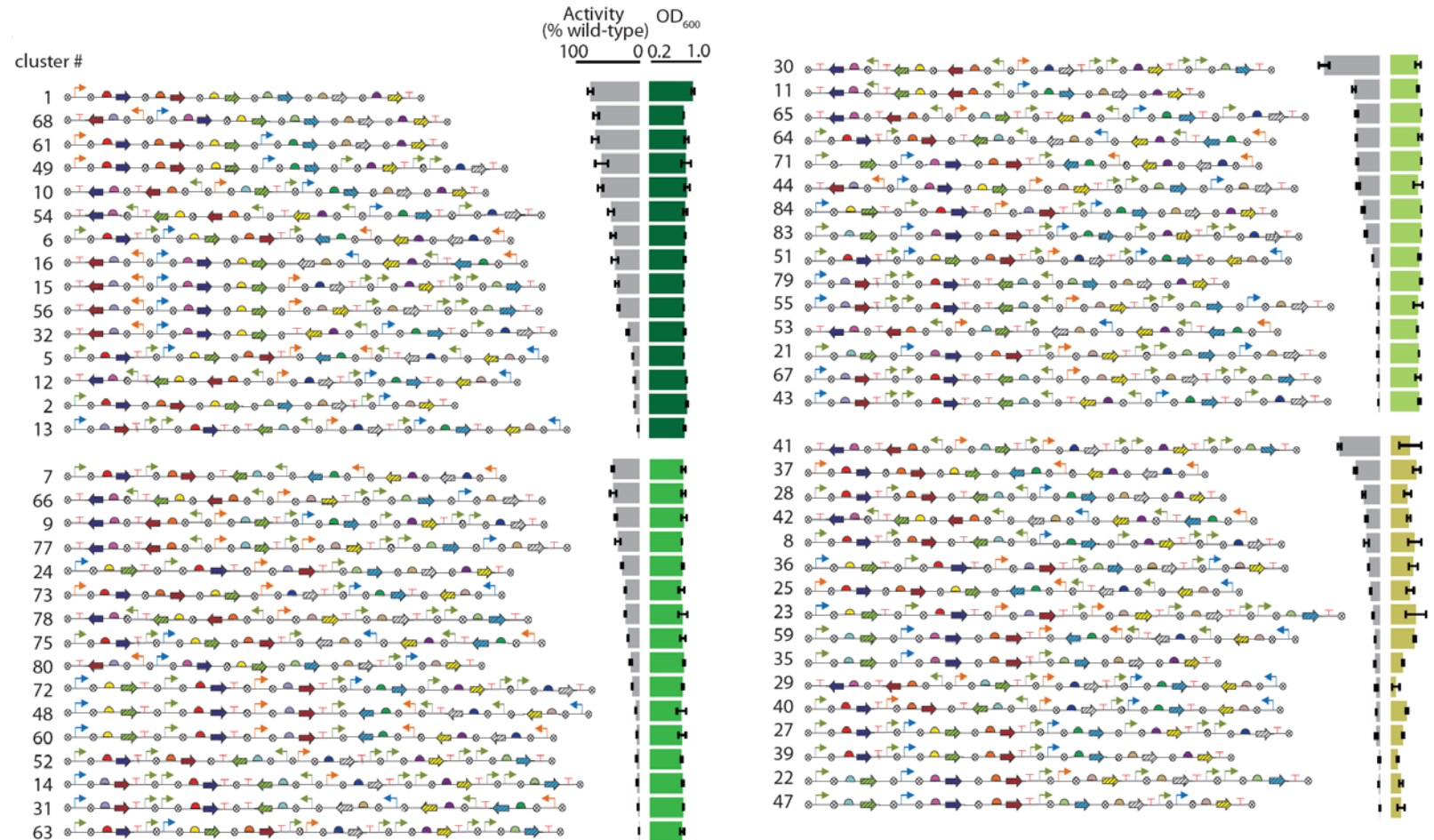
- Gene expression changes upon host transfer
- Even conservative host changes can break a system



Robert G. Egbert, and Eric Klavins PNAS 2012;109:16817-16822

# Gene Expression is Finicky but Important for Heterologous Production

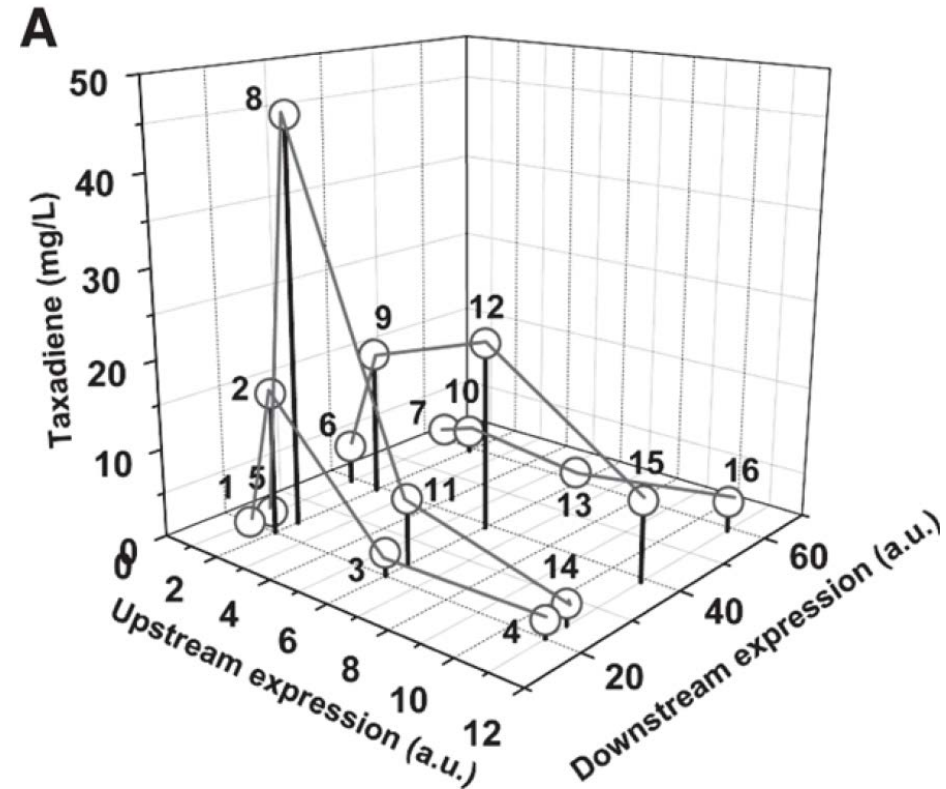
- Gene expression changes upon host transfer
- Even conservative host changes can break a system
- Permuting expression over multiple genes dramatically impacts system performance



Smanski et al. Voigt (2014) Nat. Biotechnol. 32:1241-1249

# Gene Expression is Finicky but Important for Heterologous Production

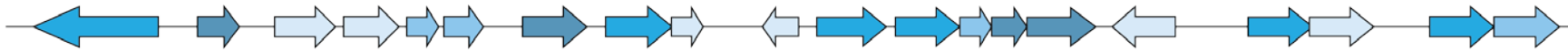
- Gene expression changes upon host transfer
- Even conservative host changes can break a system
- Permuting expression over multiple genes dramatically impacts system performance
- Optimal gene expression is nuanced and non-obvious



Ajikumar PK and Stephanopoulos G (2010) Science 330:70-74

# Approach: Reconstructing Biosynthetic Gene Clusters from Parts

Natural gene cluster



1. Remove non-coding DNA
2. Eliminate non-essential genes
3. Remove transcription factors
4. Re-design CDSs



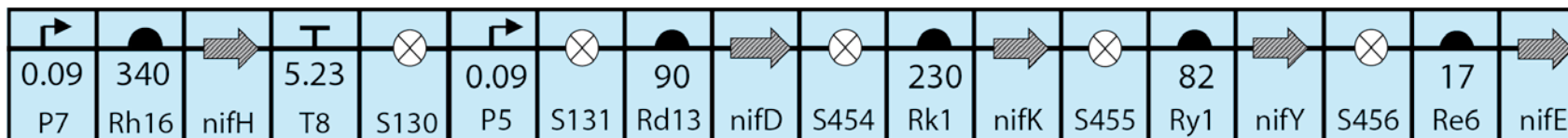
5. Clone/Synthesize genes
6. Add synthetic regulation
7. Organize into operons
8. Control with synthetic circuits



## Advantages:

- Role of every element is known
- Independent control of gene expression levels
- Ability to build and test many variant designs

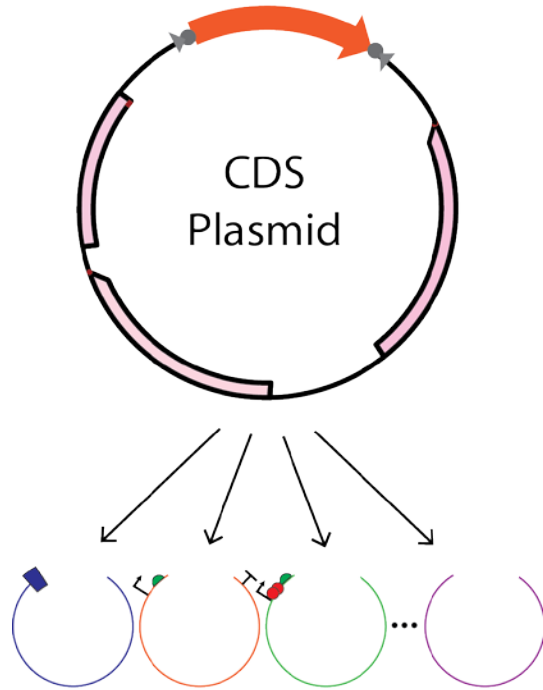
Refactored gene cluster



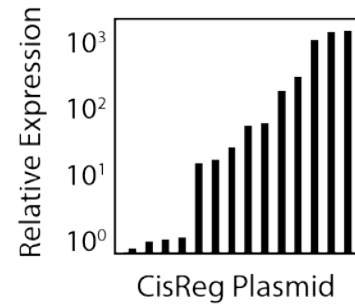
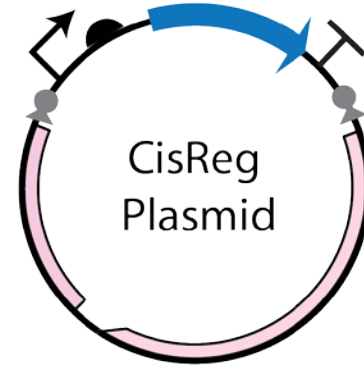


# Algorithmic DNA Assembly Pipeline: Key Features

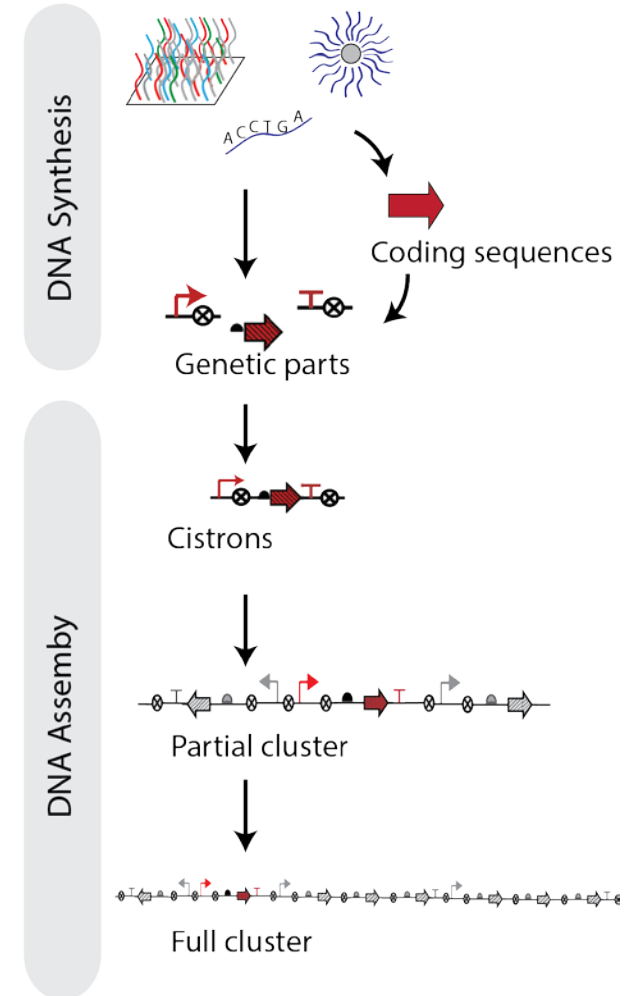
## 1. Standardized CDS Domestication



## 2. Combinatorial Quantification of Cis Regulatory Elements



## 3. Hierarchical Assembly



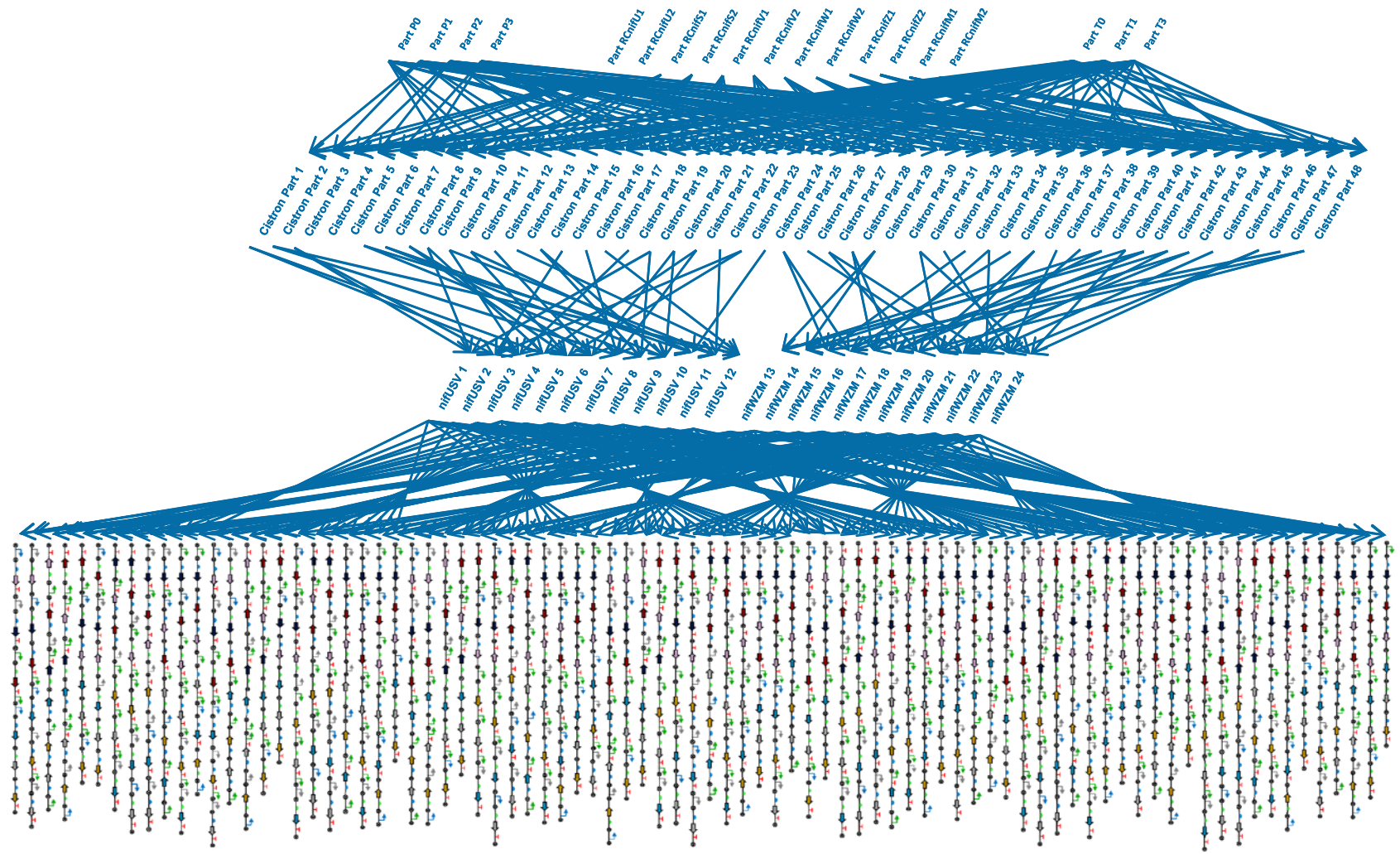
# Algorithmic DNA Assembly Pipeline For High-Throughput Plasmid Construction

Genetic Parts:

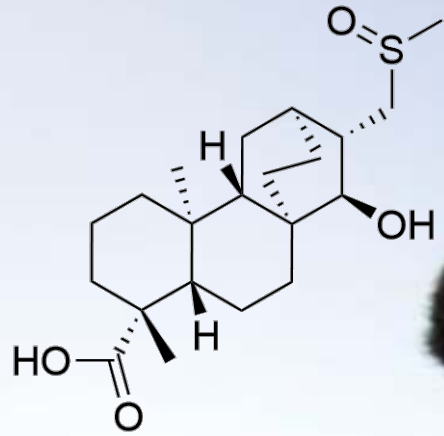
Cistron-level Constructs:

Partial Gene Clusters:

Full Gene Clusters:



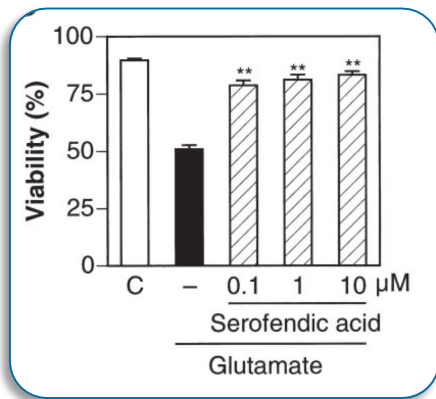
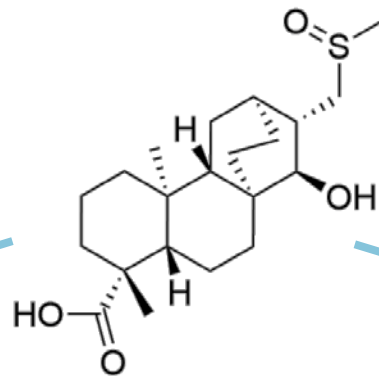
# Proof-of-concept: Biosynthesis of a Natural Product of Unknown Origin



Serofendic Acid

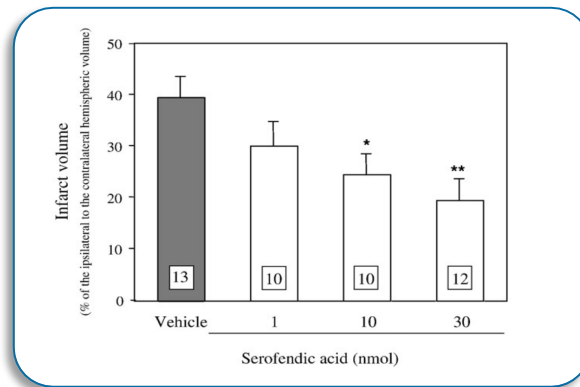


# Serofendic Acid has Diverse Therapeutic Potential



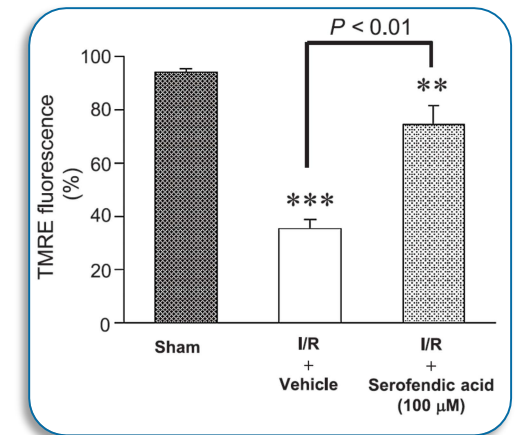
Prevents glutamate neurotoxicity

Kume, T et al. (2002) *PNAS* **99**:3288-3293.



Decreases neurological damage from stroke

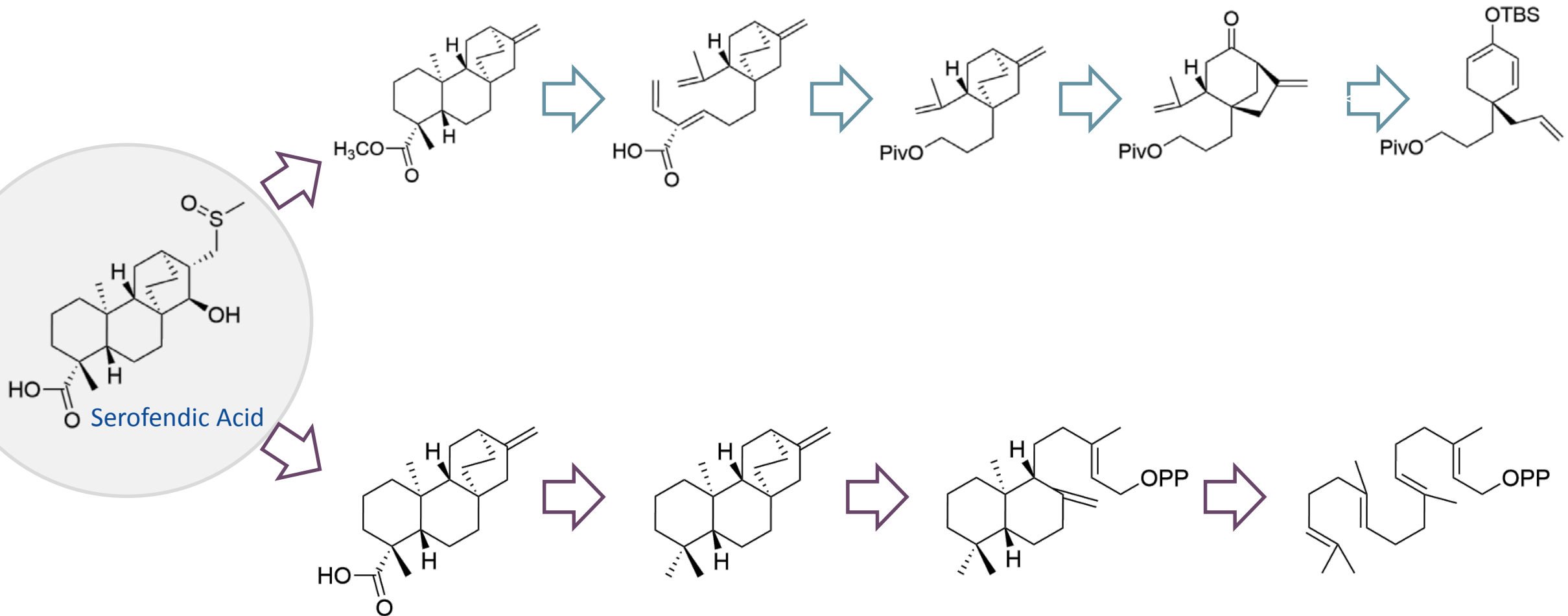
Nakamura, T et al. (2008) *Eur J Pharmacol* **586**:3288-3293.



Decreases damage from myocardial ischemia

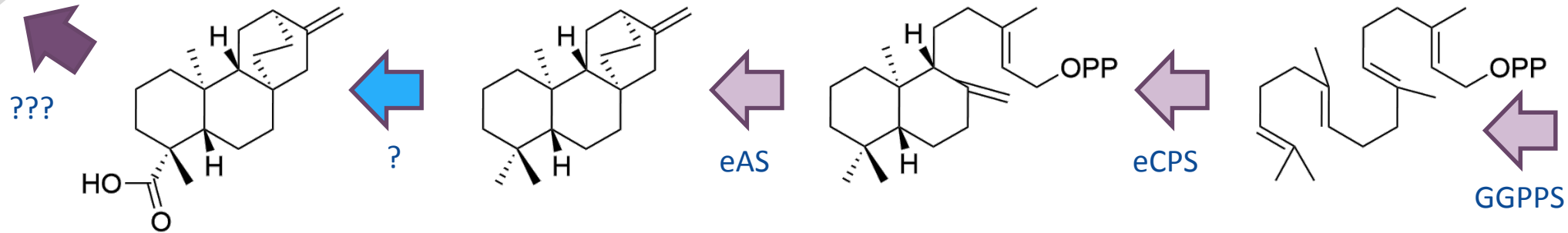
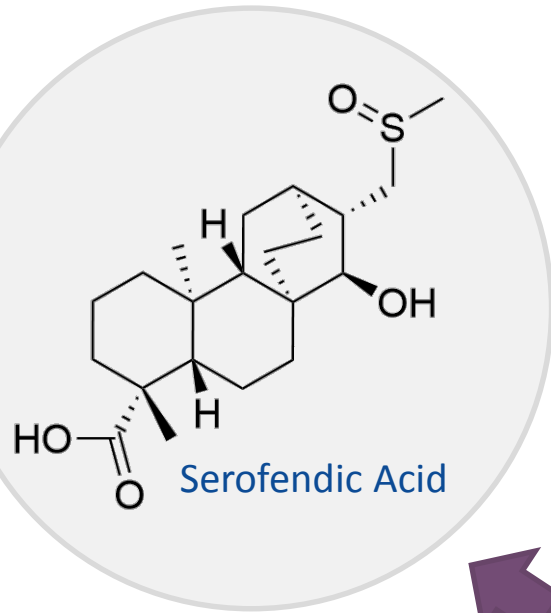
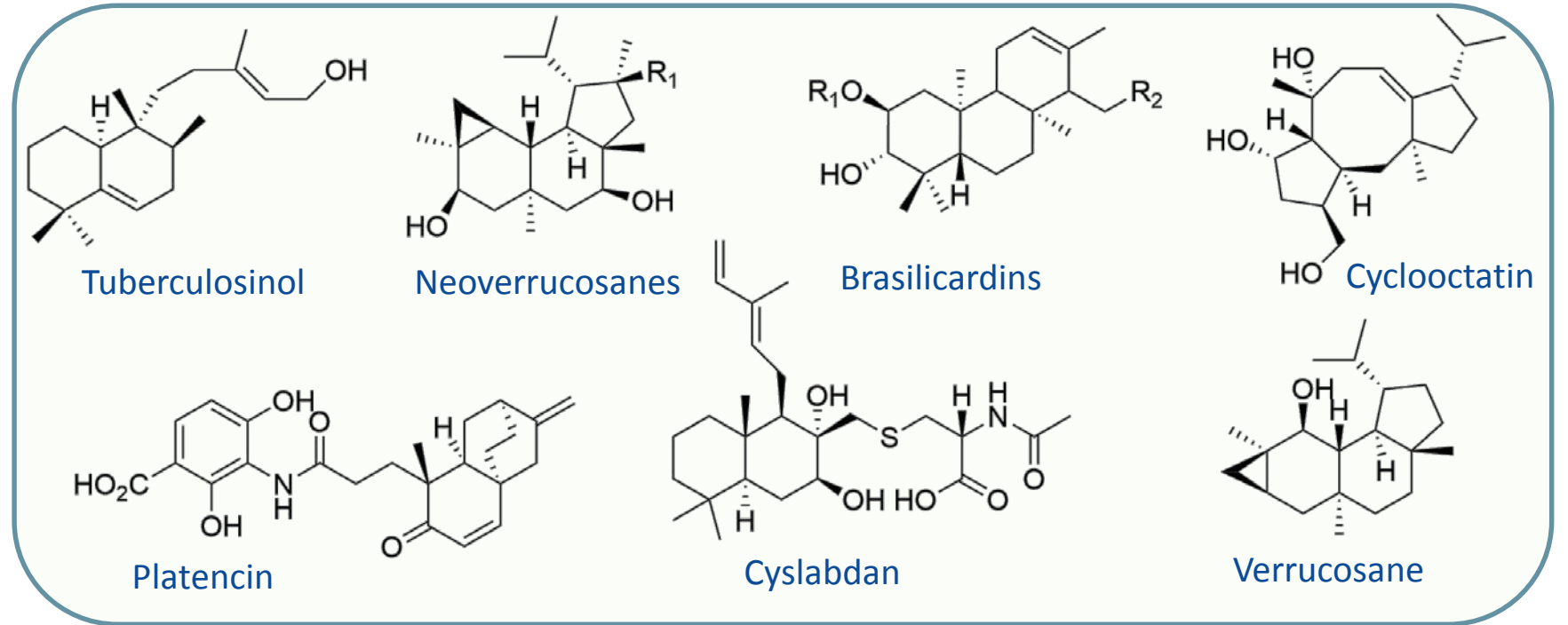
Iroji, T et al. (2014) *J Pharmacol Sci* **126**:274-280.

# Retro-Biosynthetic Approach to Design a Pathway Towards Serofendic Acid

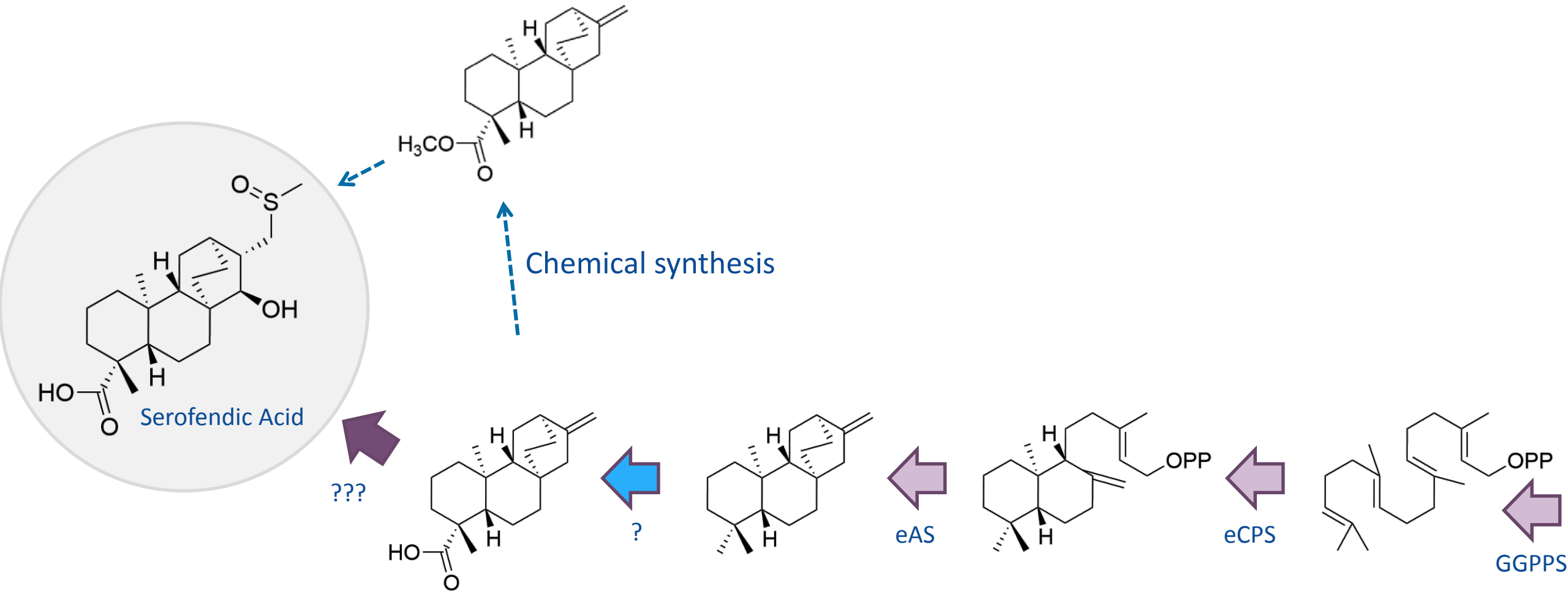


Toyota, M et al. (2005) *Org Lett* 7:3929-3932.

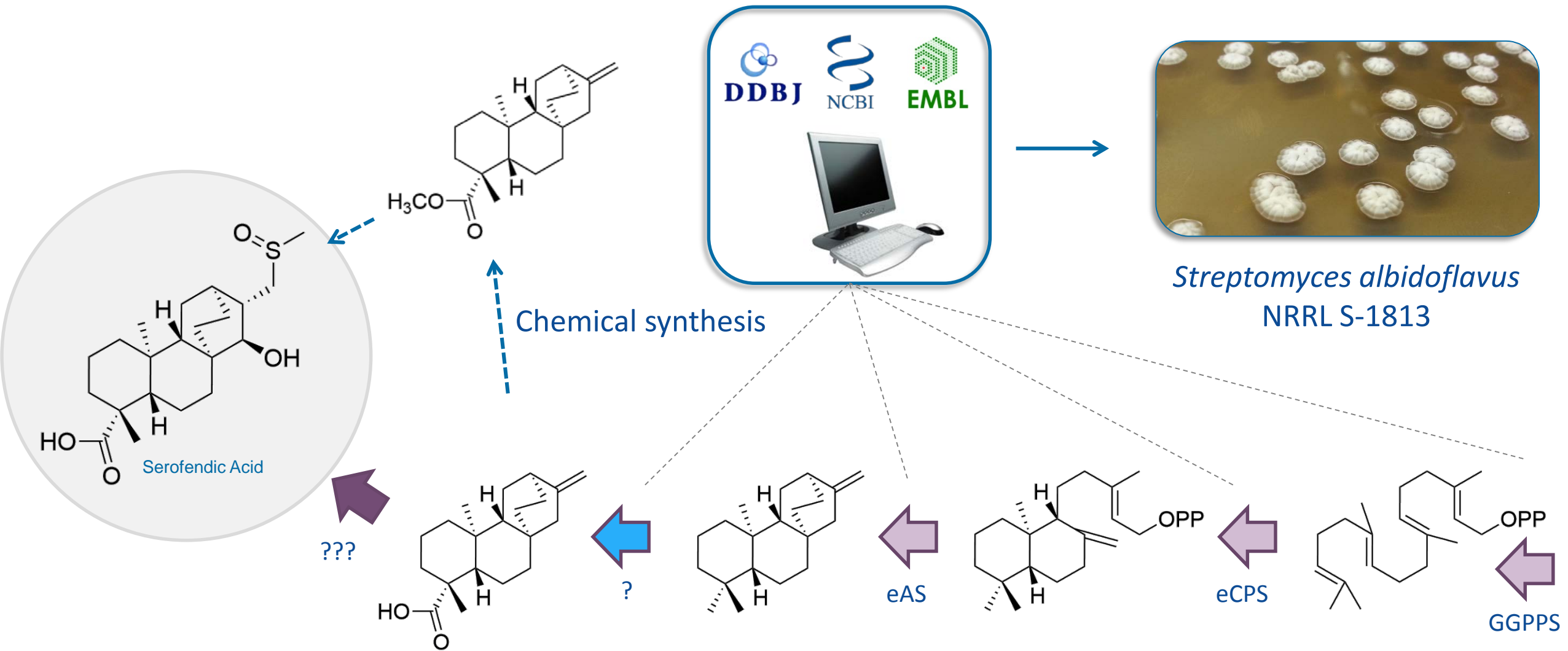
# Retro- Biosynthetic Approach to Design a Pathway Towards Serofendic Acid



# Retro- Biosynthetic Approach to Design a Pathway Towards Serofendic Acid



# Retro- Biosynthetic Approach to Design a Pathway Towards Serofendic Acid

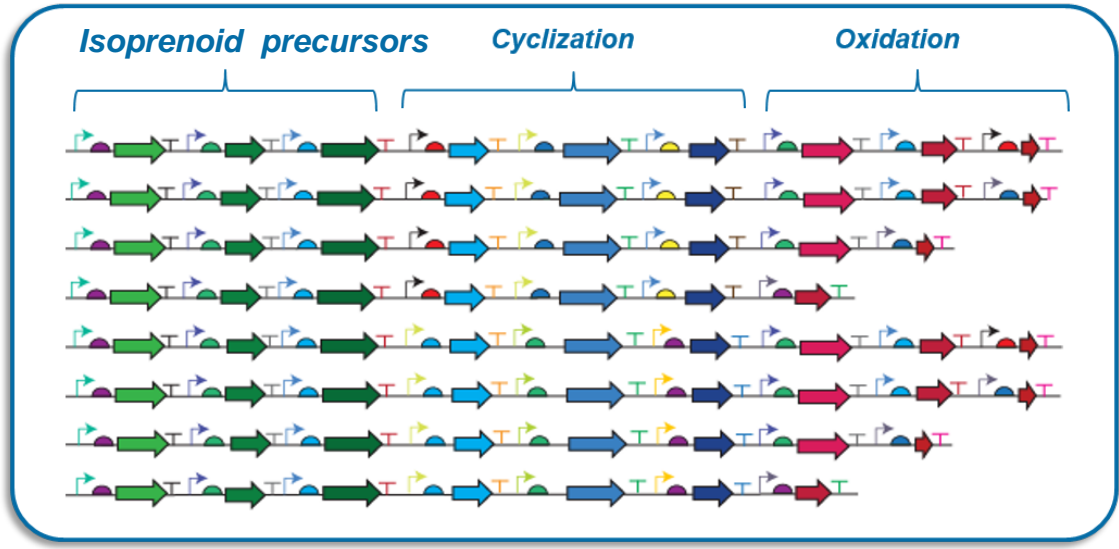
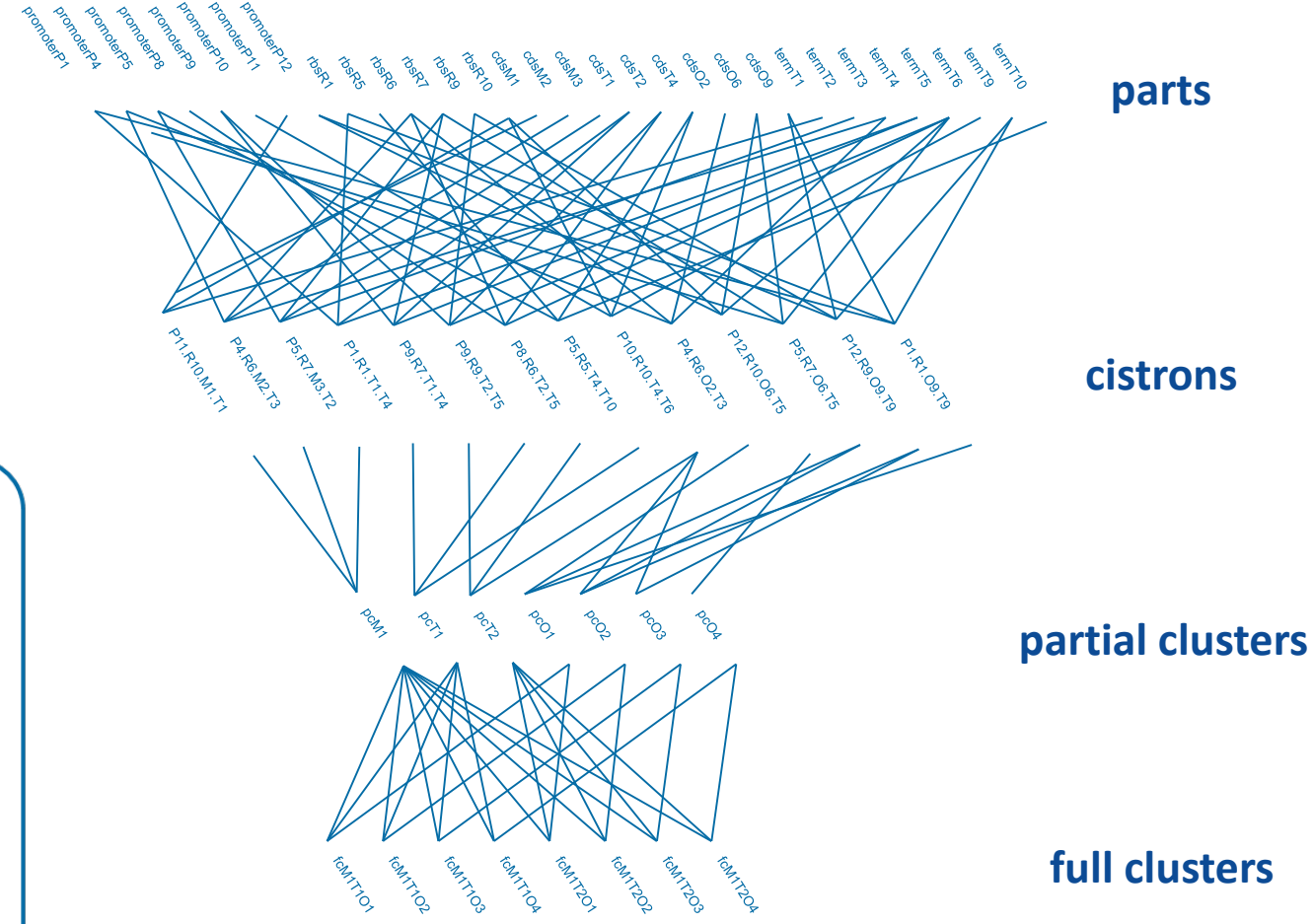
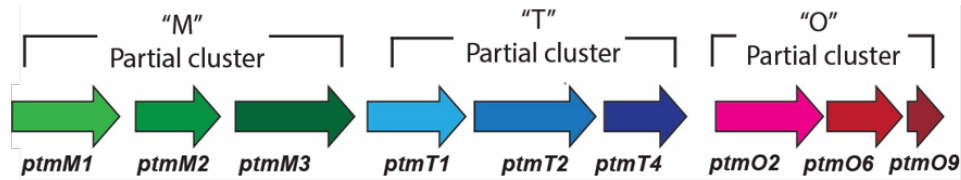




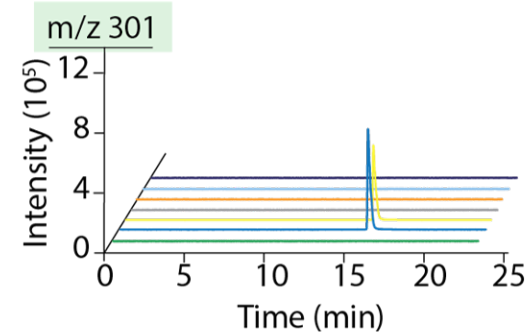
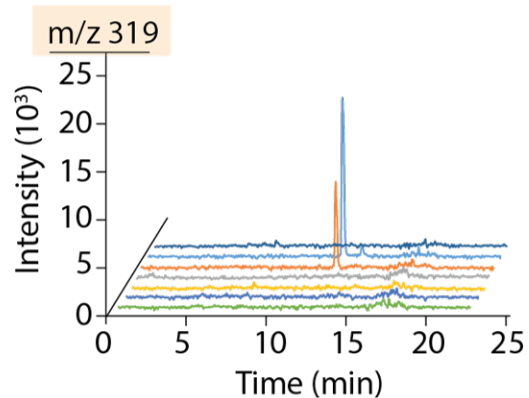
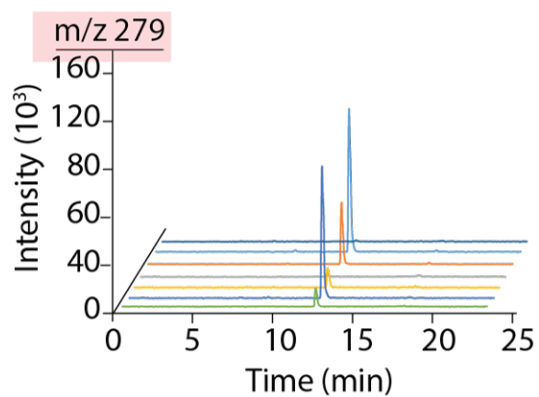
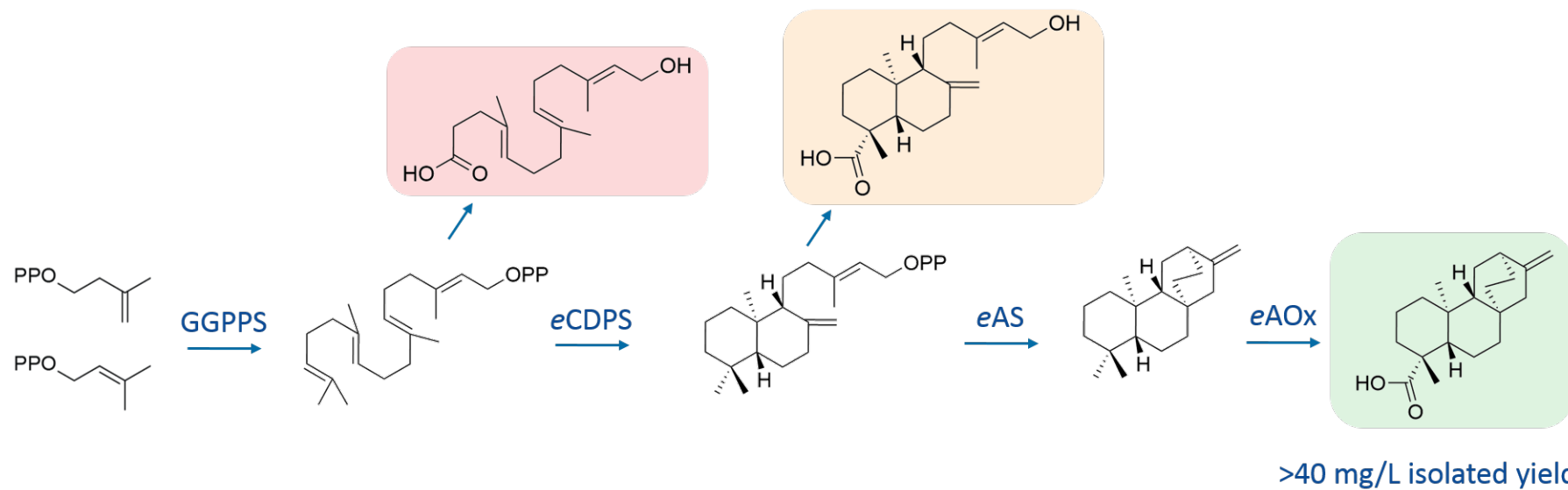
# Mini-Library Designed and Built to Screen for *Ent*-atiserenoic Acid Production

8 designs vary according to:

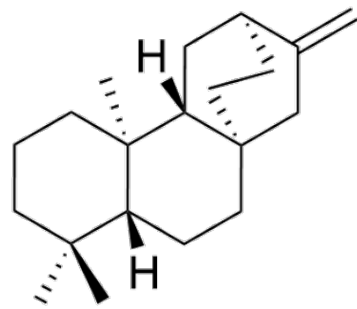
- Promoter strength
- Ribosome binding site strength
- Gene content



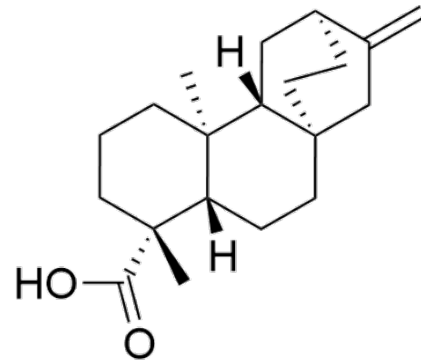
# Initial Library of Synthetic Gene Clusters Produces *Ent*-atiserenoic Acid & Congeners



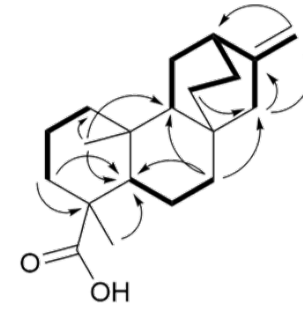
# (A) 6-Electron Oxidation of Methyl Group by P450 Monooxygenase PtnO2



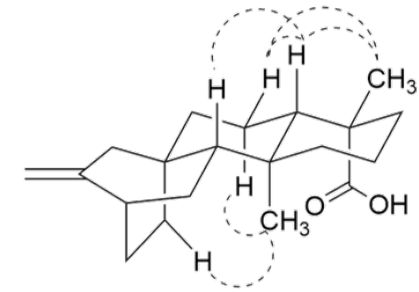
*ent*-Atiserene  
[M-H] = 271



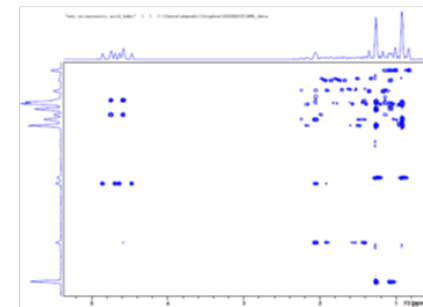
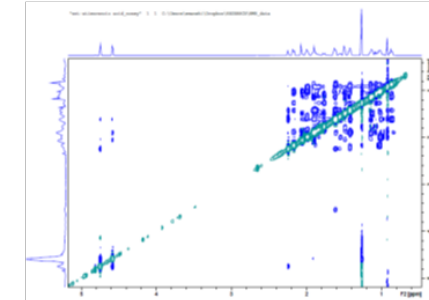
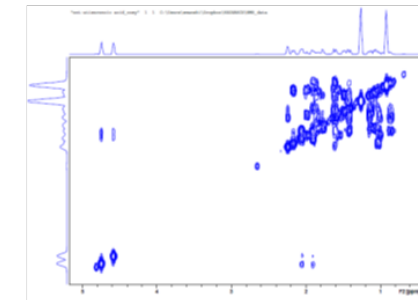
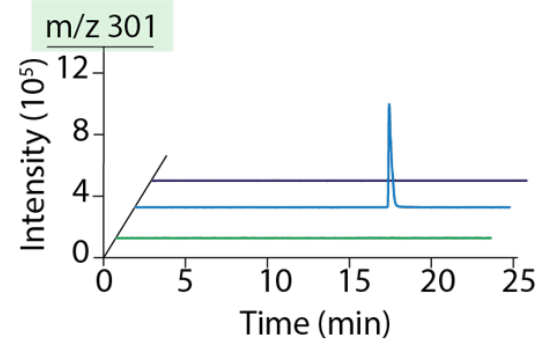
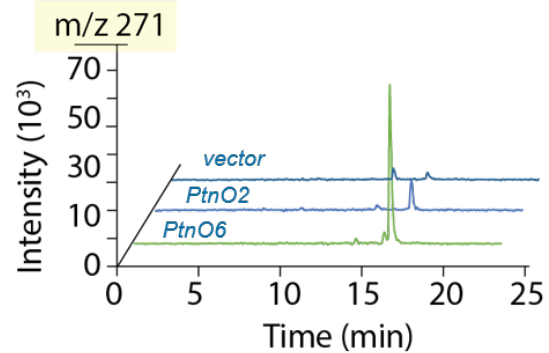
*ent*-Atiserenoic Acid  
[M-H] = 301



[COSY, HMBC]



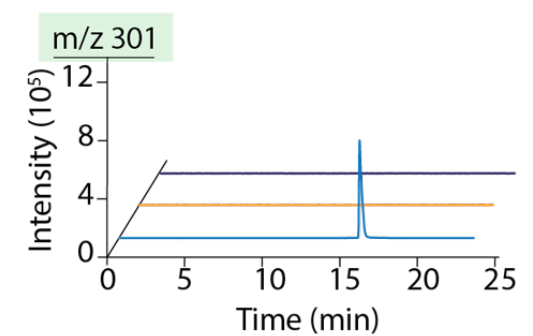
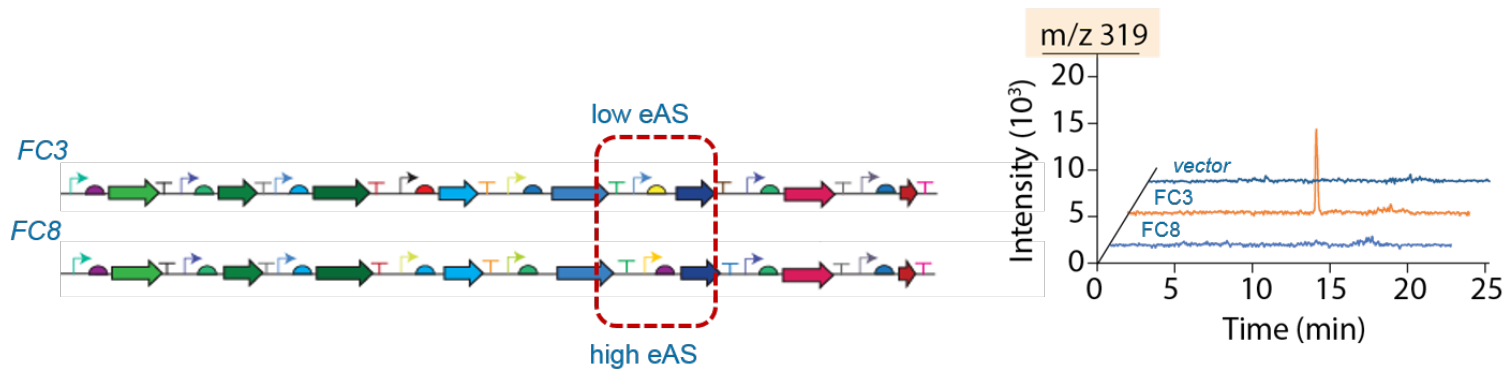
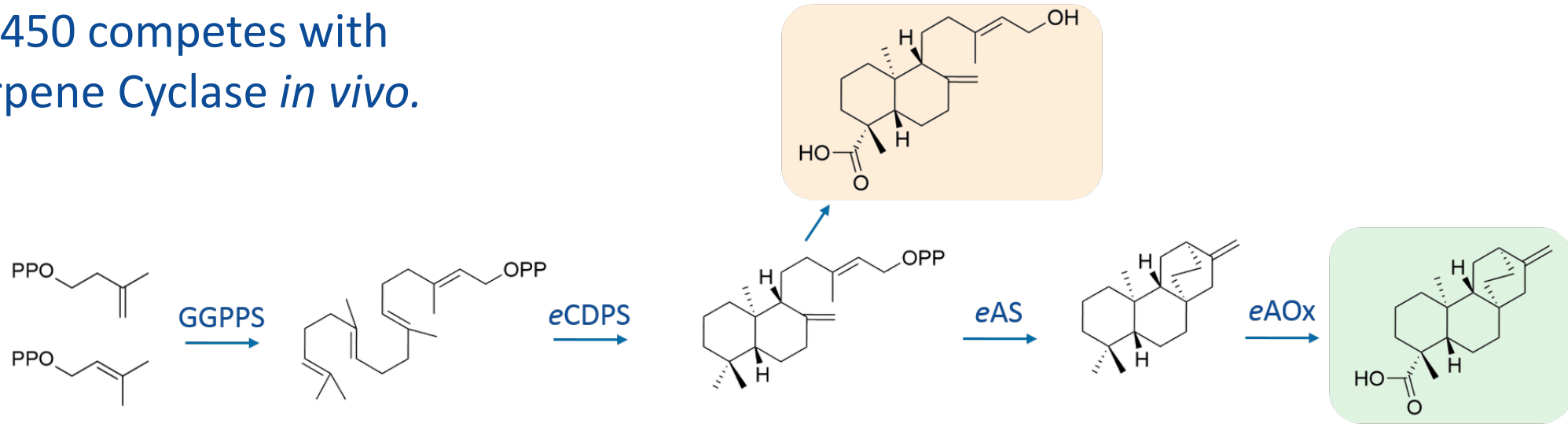
[NOESY]



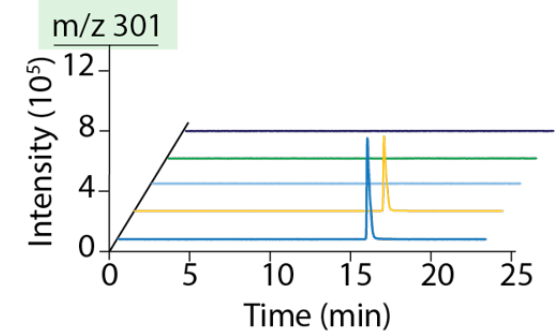
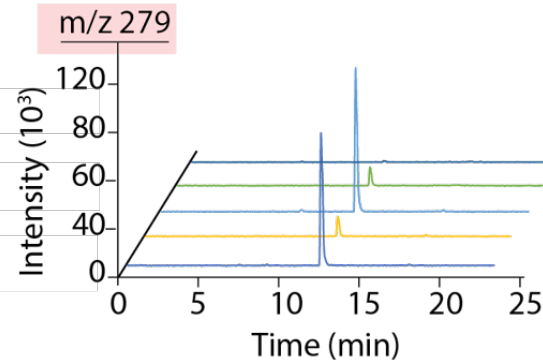
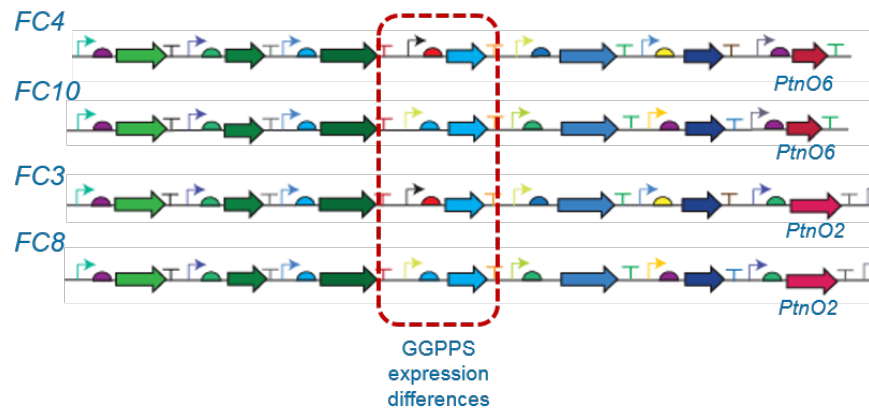
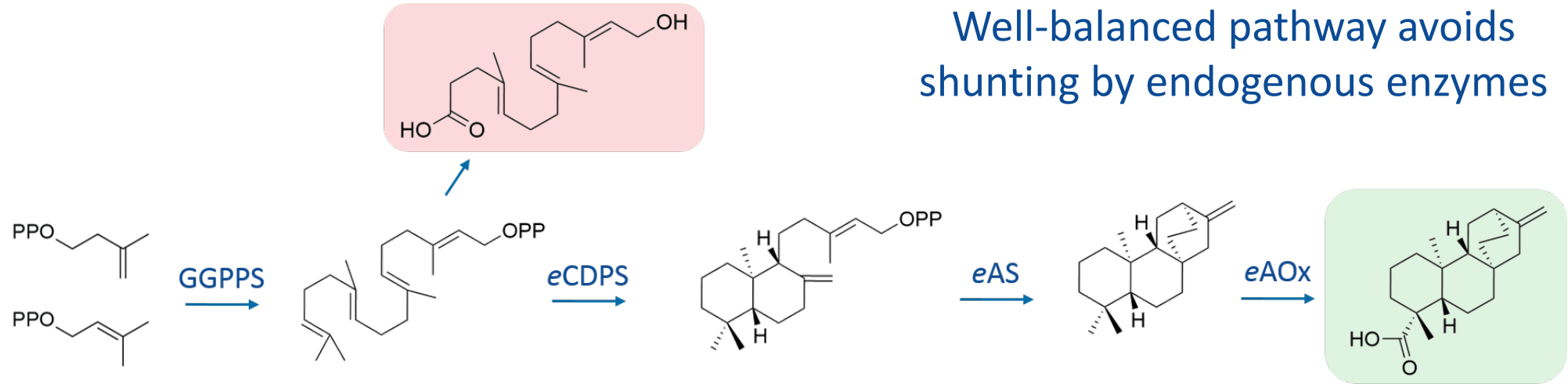
Oxidation by PtnO2, a new cytochrome P450, shows high regioselectivity.

# (B) Poorly-Tuned Gene Expression Leads to Shunt Metabolite Production

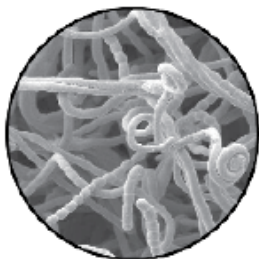
P450 competes with Terpene Cyclase *in vivo*.



# (C) Poorly-Tuned Gene Expression Allows Interference by Non-Pathway Enzymes

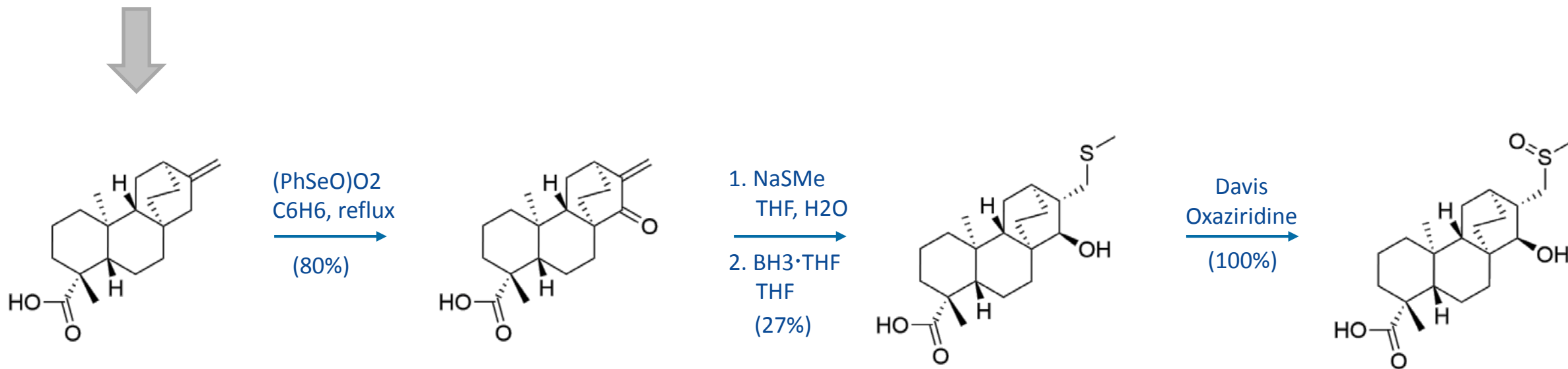


# Semi-Synthesis of Serofendic Acid and Derivatives

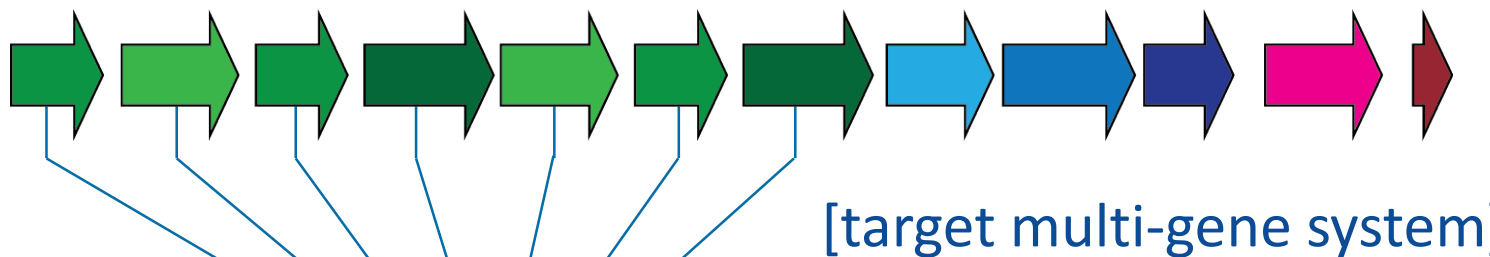


recombinant *Streptomyces*

- Initial extraction yielded 40 mg / L *ent*-Atiserenoic acid.
- Formal synthesis completed without methylester protecting group
- Semi-synthesis reduced complexity from 17 to 4 steps and increased yield ~10-fold
- Facile derivatization of *ent*-Atiserenoic acid demonstrated



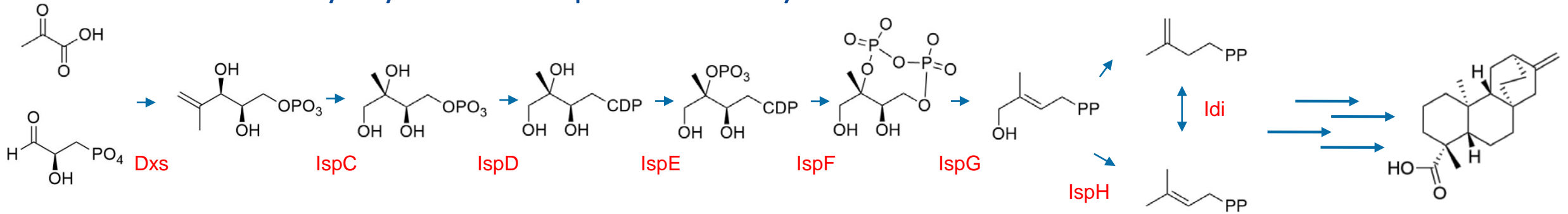
# Current Efforts are Focused on Improving Titer Through Multivariate Optimization



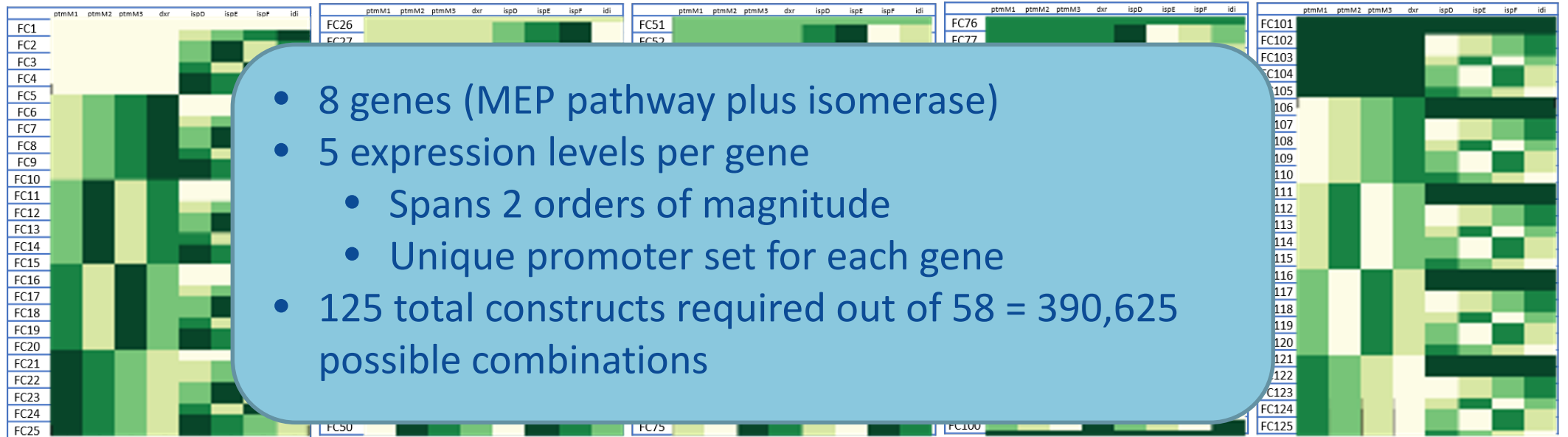
<p>Nelder-Mead simplex algorithm</p> <p>[hill climbing]</p>	<table border="1"> <tbody> <tr><td>1</td><td>+</td><td>+</td><td>+</td><td>-</td><td>+</td><td>-</td><td>-</td></tr> <tr><td>2</td><td>-</td><td>+</td><td>+</td><td>+</td><td>-</td><td>+</td><td>-</td></tr> <tr><td>3</td><td>-</td><td>-</td><td>+</td><td>+</td><td>+</td><td>-</td><td>+</td></tr> <tr><td>4</td><td>+</td><td>-</td><td>-</td><td>+</td><td>+</td><td>+</td><td>-</td></tr> <tr><td>5</td><td>-</td><td>+</td><td>-</td><td>-</td><td>+</td><td>+</td><td>+</td></tr> <tr><td>6</td><td>+</td><td>-</td><td>+</td><td>-</td><td>-</td><td>+</td><td>+</td></tr> <tr><td>7</td><td>+</td><td>+</td><td>-</td><td>+</td><td>-</td><td>-</td><td>+</td></tr> <tr><td>8</td><td>-</td><td>-</td><td>-</td><td>-</td><td>-</td><td>-</td><td>-</td></tr> </tbody> </table> <p>Plackett-Burman model</p> <p>[variable prioritization]</p>	1	+	+	+	-	+	-	-	2	-	+	+	+	-	+	-	3	-	-	+	+	+	-	+	4	+	-	-	+	+	+	-	5	-	+	-	-	+	+	+	6	+	-	+	-	-	+	+	7	+	+	-	+	-	-	+	8	-	-	-	-	-	-	-	<p>Variable 2</p> <p>Variable 1</p> <p>Central Composite design</p> <p>[interaction estimation]</p>	<p>Variable 3</p> <p>Variable 2</p> <p>Variable 1</p> <p>Full Factorial design</p> <p>[full landscape visualization]</p>
1	+	+	+	-	+	-	-																																																												
2	-	+	+	+	-	+	-																																																												
3	-	-	+	+	+	-	+																																																												
4	+	-	-	+	+	+	-																																																												
5	-	+	-	-	+	+	+																																																												
6	+	-	+	-	-	+	+																																																												
7	+	+	-	+	-	-	+																																																												
8	-	-	-	-	-	-	-																																																												

# Current efforts are focused on improving titer through multivariate optimization

## Methylerythritol Phosphate Pathway

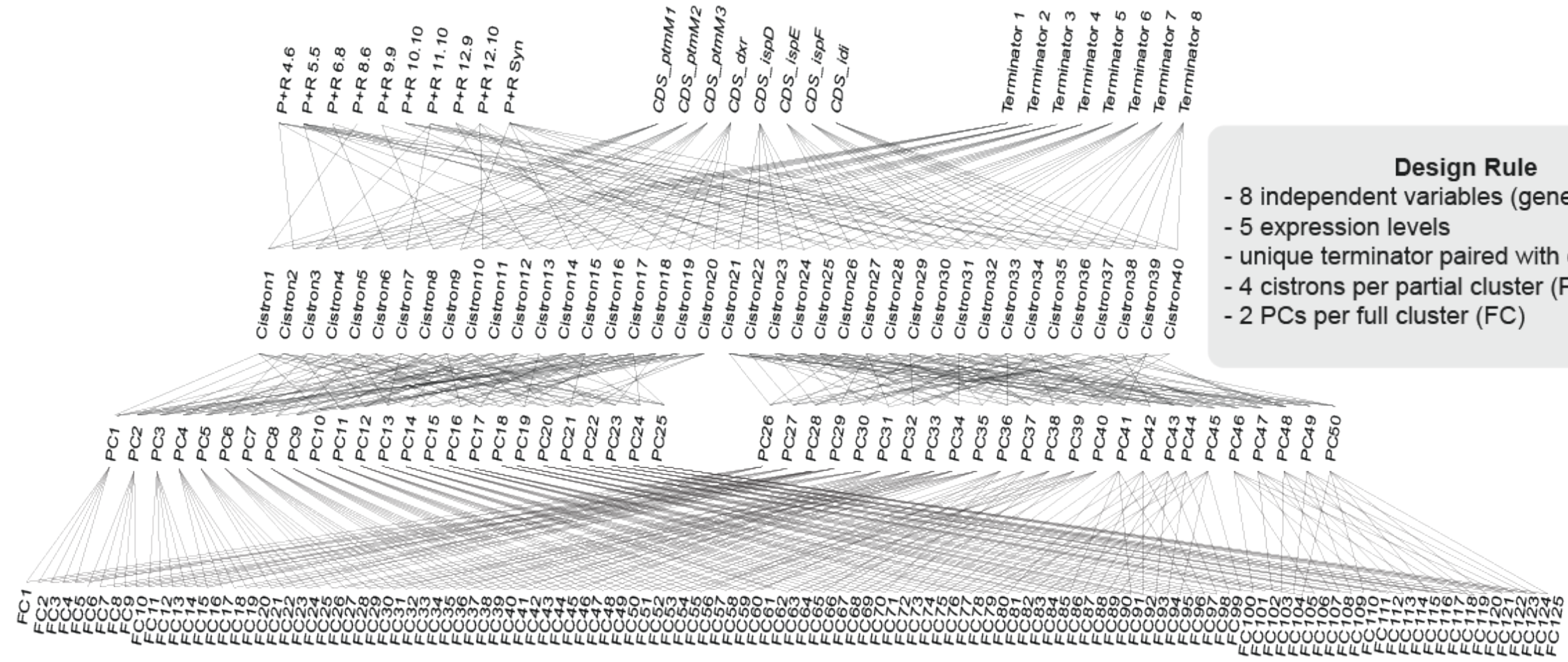


## 5-level fractional factorial design:





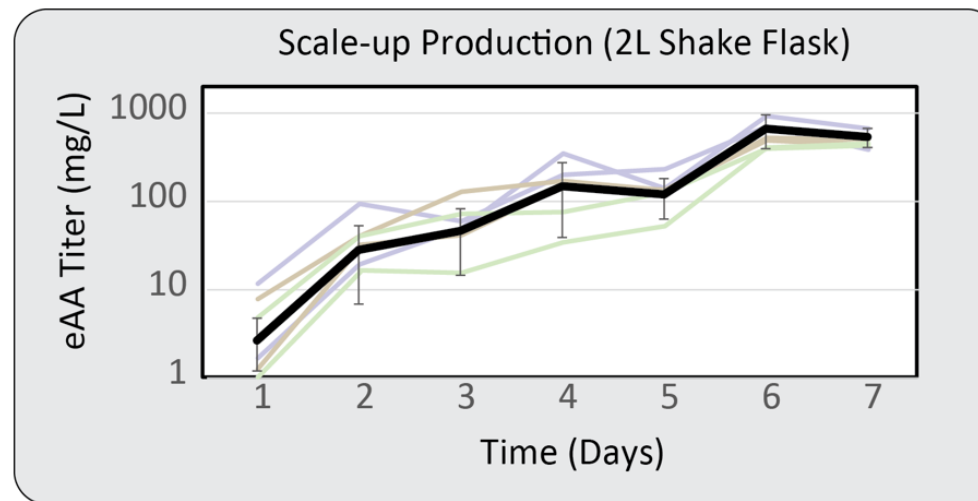
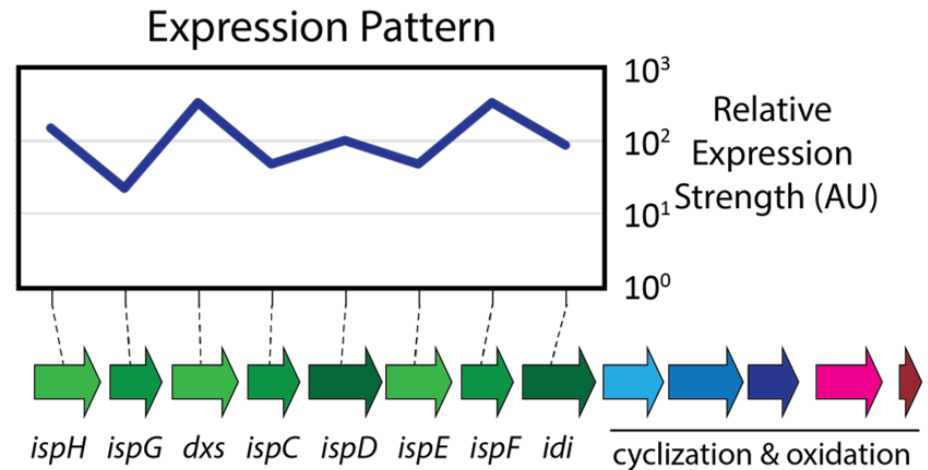
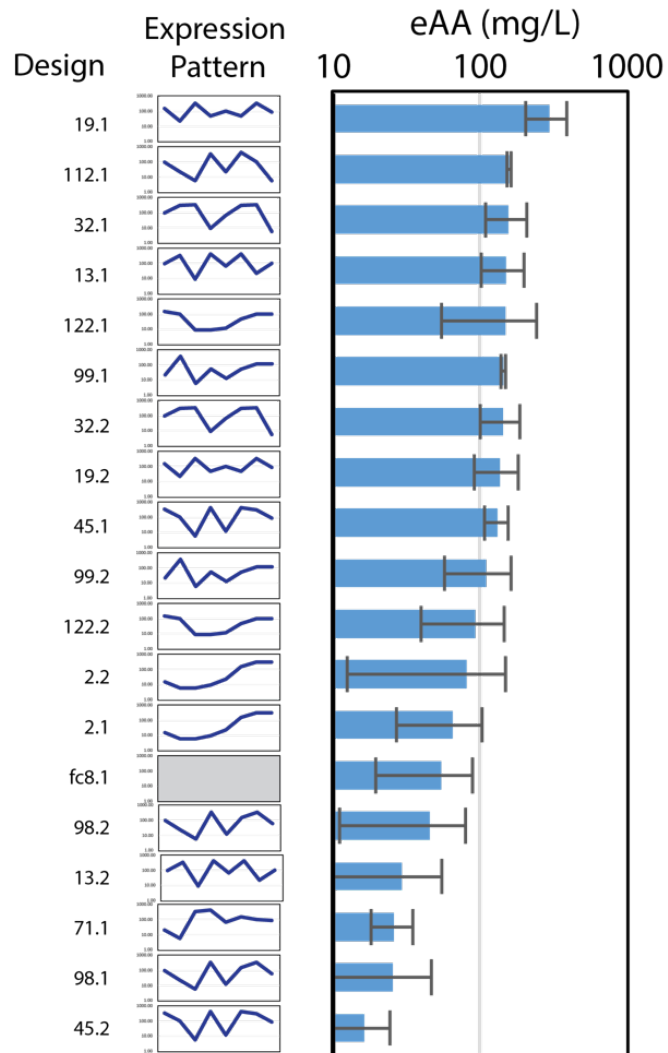
# Current Efforts are Focused on Improving Titer Through Multivariate Optimization



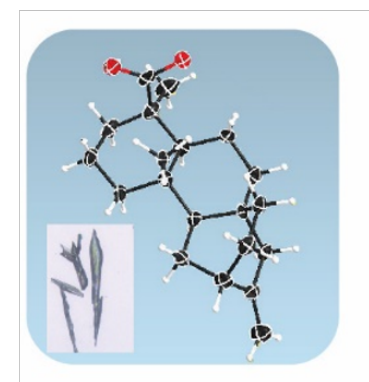
**Design Rule**

- 8 independent variables (gene expression)
- 5 expression levels
- unique terminator paired with each rCDS
- 4 cistrons per partial cluster (PC)
- 2 PCs per full cluster (FC)

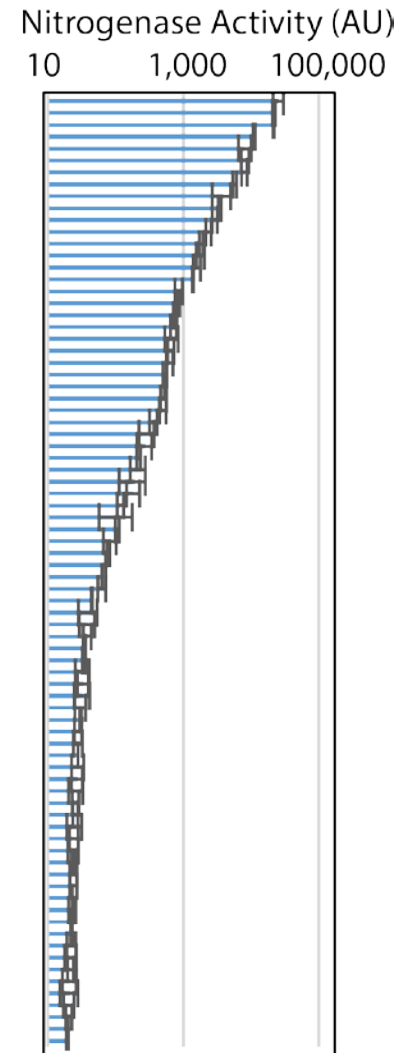
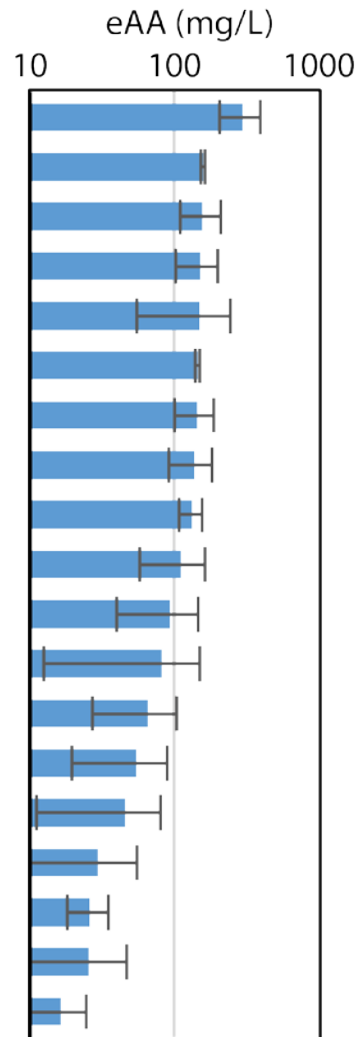
# Preliminary Results Show Several Improved Strains



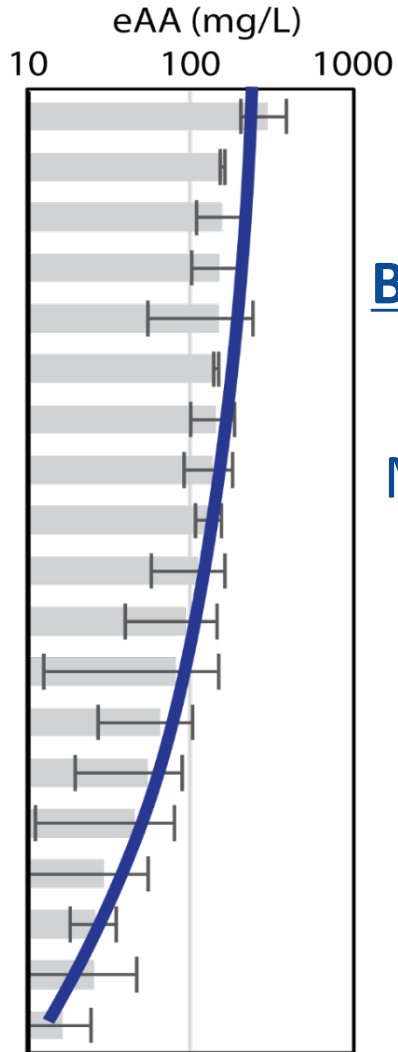
Recent titers  
>500 mg/L



# Preliminary Results Show Several Improved Strains



# Robustness to Varying Expression Levels Will Not Be the Same for Every System



## Diterpene Biosynthesis

Secondary Metabolism

Discrete Enzymes

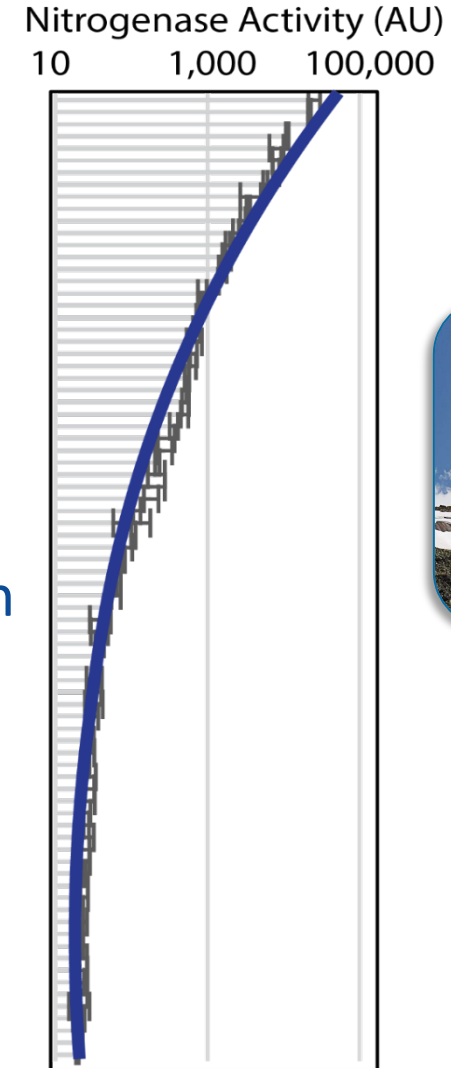
vs.

## Nitrogen Fixation

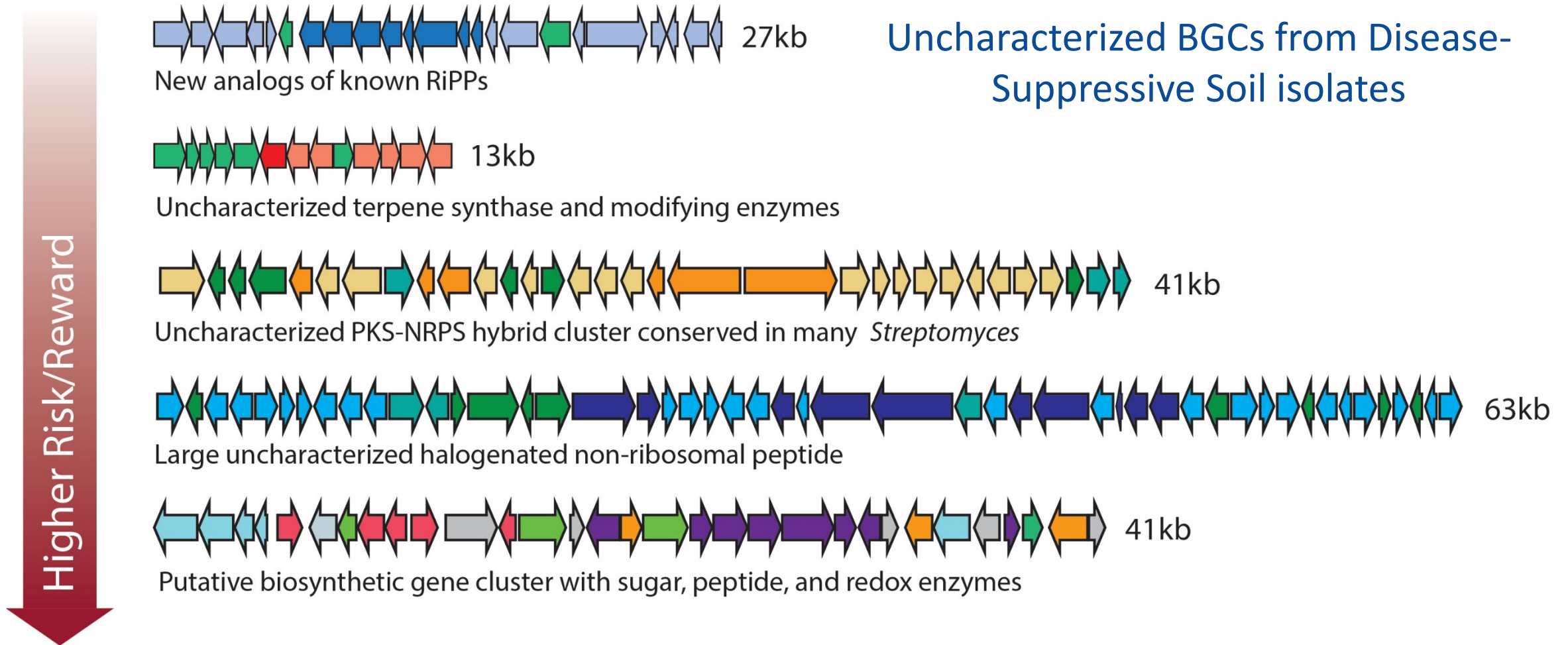
Primary Metabolism

Multi-protein Complexes

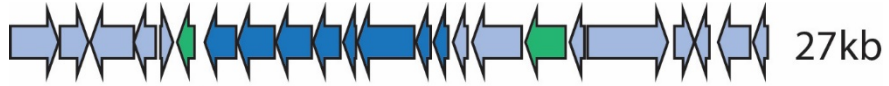
vs.



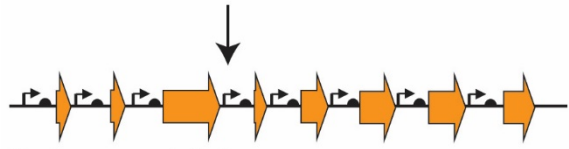
# Extending the DNA Assembly Pipeline for New Compound Discovery



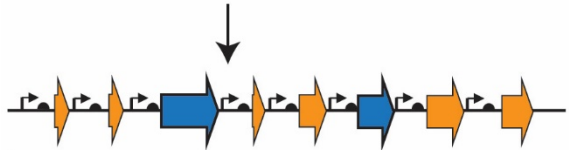
# Family of Putative Thiazole-Containing Molecules



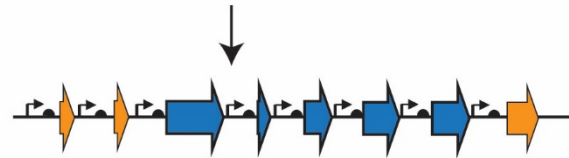
New analogs of known RiPPs



Refactored 1.0



Refactored 2.0

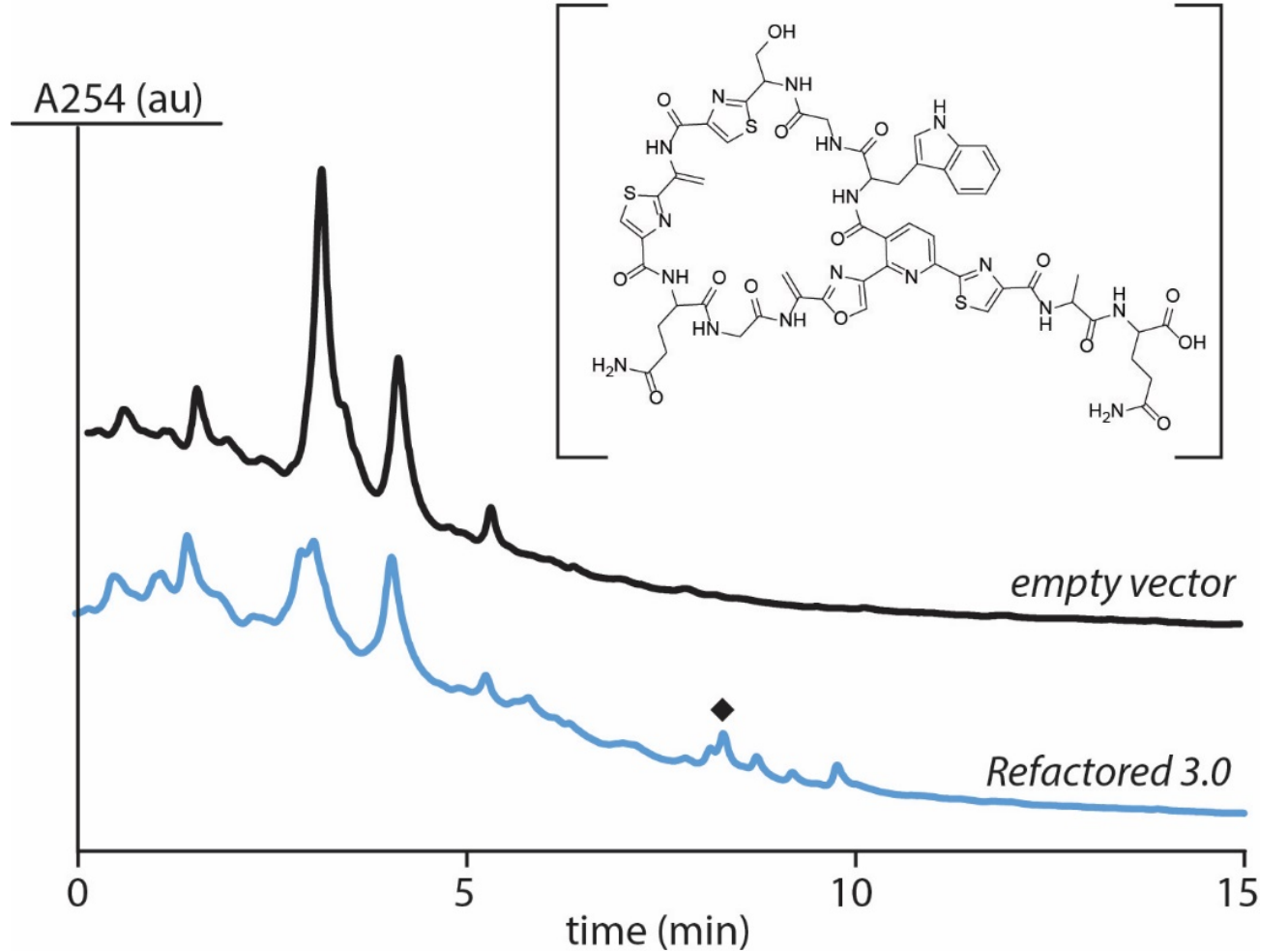


Refactored 3.0

Wildtype CDS

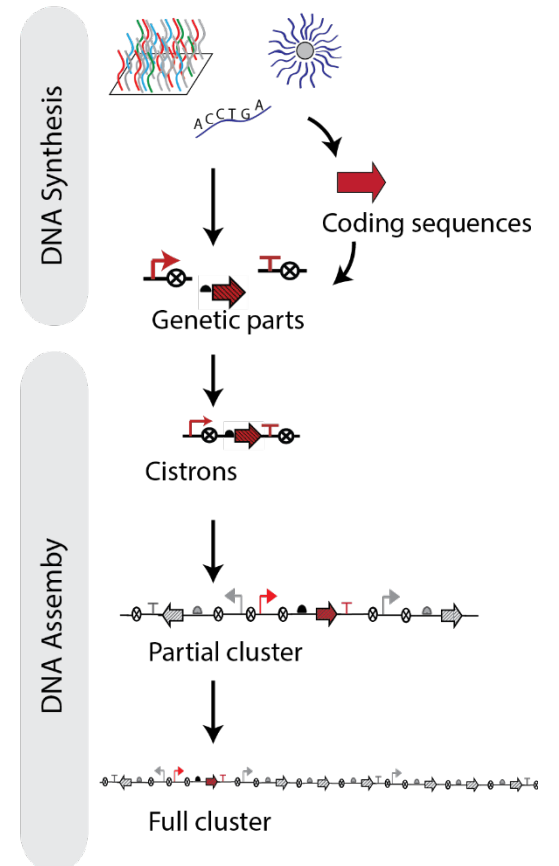
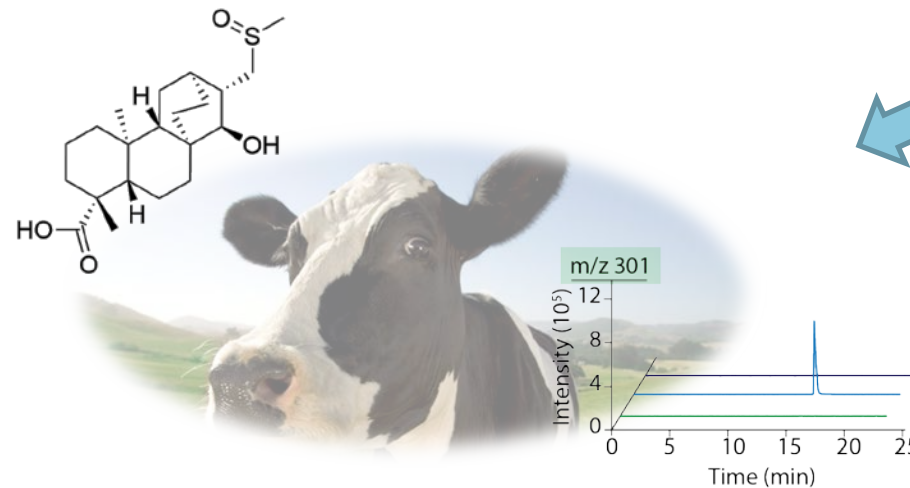
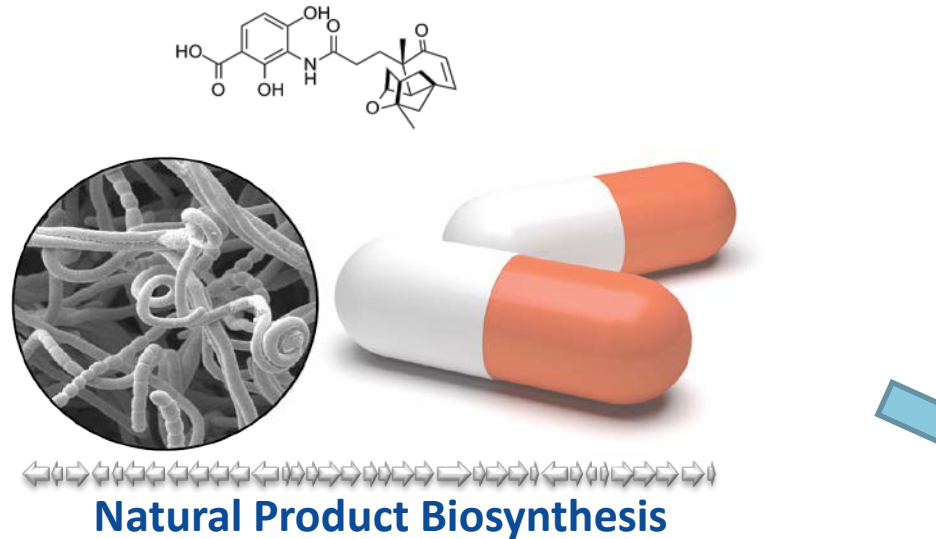
Codon altered CDS

Thiazole BGC with proposed role  
in interspecies signaling.



# Summary: DNA Assembly for Engineering Natural Product Biosynthesis

- Independent control of gene expression is important for high-titer heterologous production
- Developed an algorithmic DNA assembly pipeline compatible with *Streptomyces*
- Demonstrated a new sustainable route to Serofendic acid, a diterpenoid natural product of unknown origin



## Collaborators

**Linda Kinkel (UMN)**

Dr. Lindsey Hansen

Dr. Zewei Song

Dan Schlatter

## Funding and Support

BTI Biocatalysis Grant

UMN Futures Grant

Damon Runyon Cancer

Research Foundation

Joint Genome Institute

DARPA



Dr. Maciej Maseko  
Dr. Christopher Stach

**Dr. Dimitri Perusse**

**Suzie Hsu**

Stephen Heinsch

Carolyn Malecha

Matt Zinselmeier

**Thomas Hougard**

Mariela Rivera-DeJesus

Blake Everett



# GenScript: Your Reliable Research Partner

- Serving life scientists for **more than 15 years** as a **leading CRO** (Contract Research Organization) offering a variety of services, reagents and products
- Headquarters in **Piscataway, NJ**
- Facilities and branches in **China, Europe & Japan**
- The **number one** gene synthesis provider in the world
- **30,000** customers in **90+ countries**
- Recipient of **CRO Leadership Award** in productivity and regulatory affairs
- **≥30,000** journal article citations
- An **iGEM** partner
- The **only commercial entity** invited to participate in the Synthetic Yeast Genome project (**Sc2.0**)
- A member of **GP-write advisory board** due expertise in gene synthesis



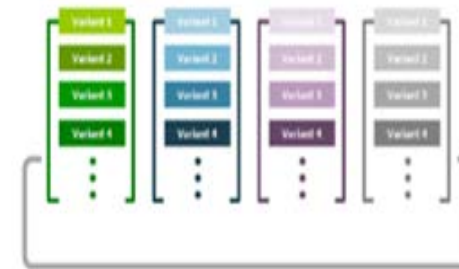
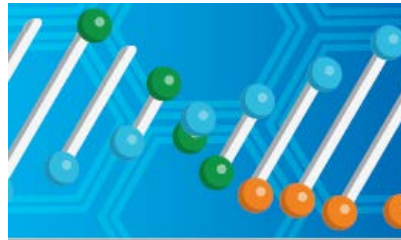
# Let DNA Building Experts Speed Up Your Metabolic Pathway and Microbial Strain Engineering Process!

## Synthetic Biology Portfolio

Gene Synthesis

DNA Fragments

Combinatorial  
DNA Libraries



Fast, Reliable & Efficient

# Partnering with Us

## Worldwide Leader

Benefiting from 15+ years of expertise in complex DNA synthesis & assembly

## Comprehensive Services

Easy access to highly-customizable & high-throughput industrial grade enabling tools

## Fast & Efficient

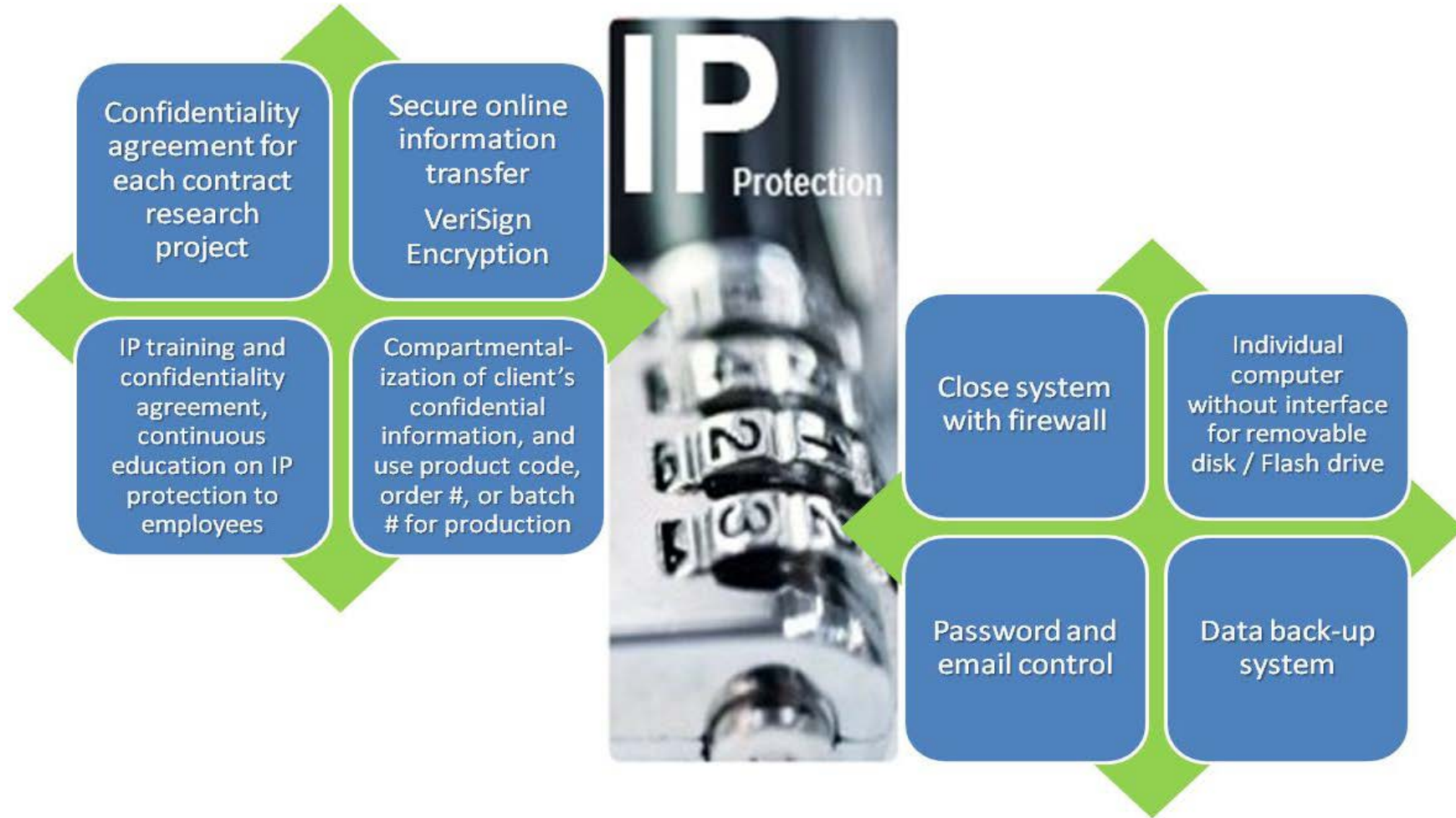
Speed up the build phase of your development cycle with time/cost-efficient options

## Trusted

Join the long list of leading life scientists in academia & industry who have advanced their synthetic biology projects with us.



# Partnering With Us: IP Protection



# Thank you!

For any questions,

please visit:

[www.genscript.com/synthetic-biology.html](http://www.genscript.com/synthetic-biology.html)

or email:

[William.tan@genscript.com](mailto:William.tan@genscript.com)  
[gene@genscript.com](mailto:gene@genscript.com)



Gene



Peptide



Protein



Antibody



Discovery



Catalog Products