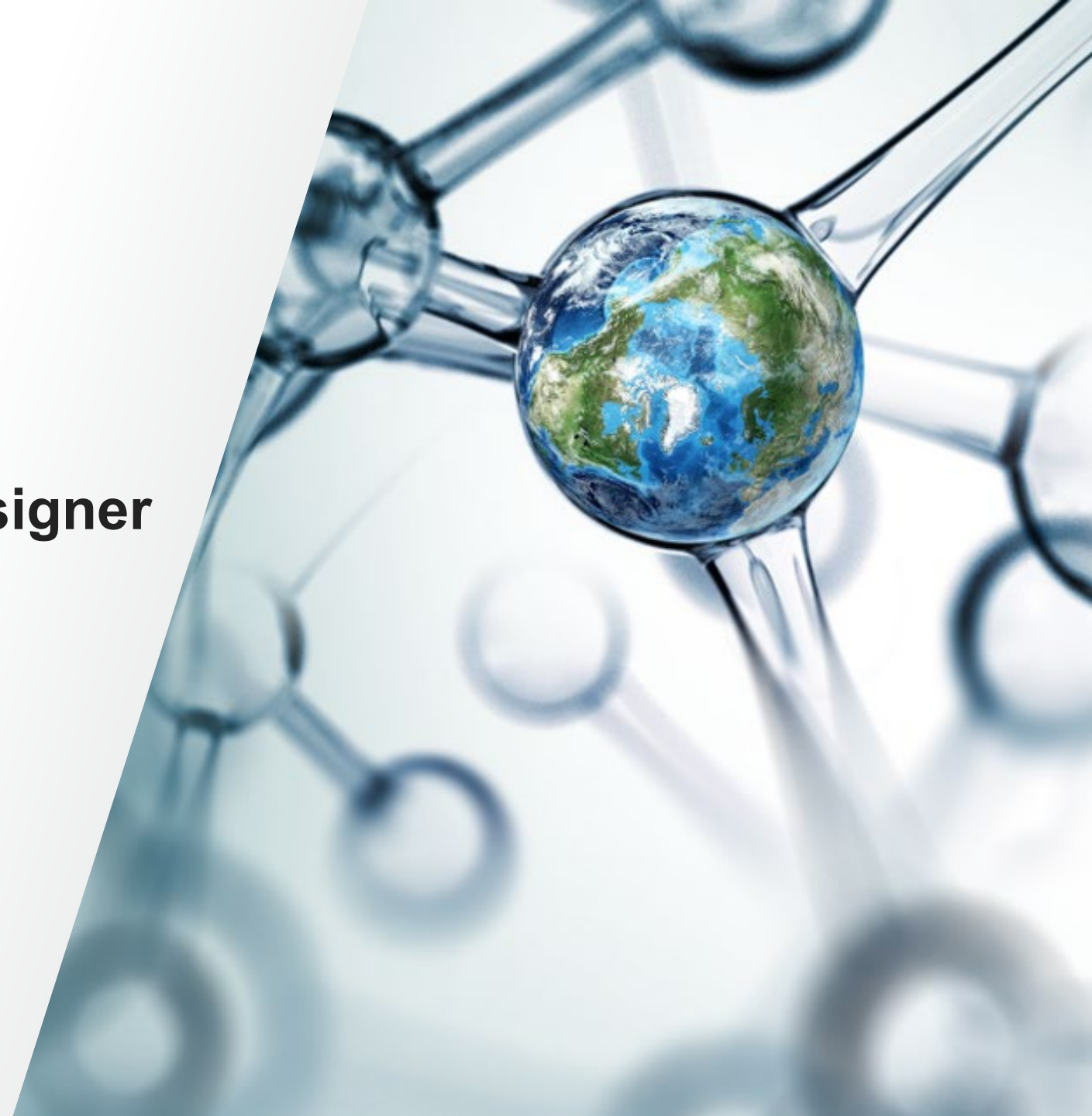


Invitrogen GeneArt Instant Designer オーダークイックガイド

バイオサイエンス事業部

 The world leader in serving science



Invitrogen™ GeneArt™ Instant Designer オーダークイックガイド

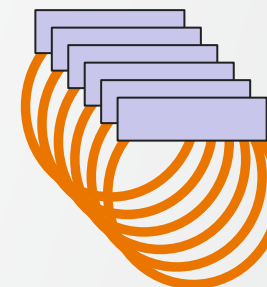
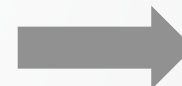
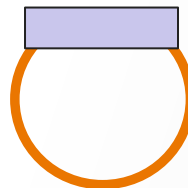
こちらは以下の内容の Web 入力クイックガイドです。(塩基配列の入力からプラスミド精製入力)

塩基配列入力

ベクター導入

プラスミド精製

ATGGTGAGC...



- 詳細なマニュアルもございますので、そちらもご参照ください。
<https://assets.thermofisher.com/TFS-Assets/BID/Reference-Materials/geneart-instant-designer-order-manual-ja.pdf>
- クイックガイドになりますので、一部の説明を省略しております。
※Webの仕様は予告なしに変更する可能性があります。

販売条件

ライフテクノロジーズ標準販売条件

<https://assets.thermofisher.com/TFS-Assets/CORP/Licensing-Information/lt-standard-sales-conditions-japan.pdf>

遺伝子合成販売条件

<https://assets.thermofisher.com/TFS-Assets/BID/Licensing-Information/gene-synthesis-condition-of-sales-en-ja.pdf>

ライフテクノロジーズ・コーポレーション研究用受託サービスに関する条件

<https://assets.thermofisher.com/TFS-Assets/BID/Licensing-Information/life-technologies-custom-work-terms-conditions-ja.pdf>

ログイン後の画面

Cloned Genes の Start New Project をクリック

What can we offer you?



Cloned Genes

(100-12000 bp)

Start New Project

[Learn more](#)



DNA Fragments

(200-3000 bp)

Start New Project

[Learn more](#)



Reorder Plasmid

Order stored plasmid

Order Now

[Learn more](#)



Custom Proteins

Expressed and purified

Request a Quote

[Learn more](#)



DNA Libraries

Directed Evolution

Request a Quote

[Learn more](#)

1. Project 名を記入(例では Example1)
2. 配列入力方法を選択(ファイルアップロード or 手入力)

1. Enter project information (Optional) Default project name will be given if not specified

Example1	Optional description
---------------------	---------------------------------

2. Select sequence input method [Download Excel template](#)



Upload Sequence File

FASTA, GenBank or Excel format



Manual Creation

Copy & paste your sequences directly
into Instant Designer

配列入力(手入力の場合)

1. 配列名を記入(例では Example1)
2. 配列を入力(コピー & ペースト可、アミノ酸配列可)
3. Continue をクリック
4. ポップアップが出たら、Continue to configure をクリック

Project: Example1



1 Sequences Delete

Back

Continue

Sequence Name	DNA or Protein Sequence Important Information for Cloning and Protein Sequence(s)	5' Region (Optional)	3' Region (Optional)
☐ Example1	☐ ATGGTGAGCAAGGGCGAGGAGCTGTTACCGGGGTGGTGGCCATCTGGTCGAGCTGGACGGCGACGTA...	Click to enter 5' region	Click to enter 3' region
☐			
☐			
☐			
☐			
☐			

ポップアップ

Everything looks good!

All uploaded sequences have passed our initial checks and are ready to import. You can continue with your configurations on individual sequences in the next step. If you would like to apply same configuration on **multiple sequences** like Optimization or Cloning parameters, you can do this by selecting the sequences and click on [Configure Selected](#).

Stay on this page (or) Continue to configure

エラー表示が出る場合や
配列追加したい場合は
引き続きこのページで操作
(配列追加は後の工程でも可能)

Review画面

1. クローニングするベクターを選択
2. 使用する制限酵素ペアを選択
(制限酵素サイトが合成配列に自動的に付加されます)
3. Edit 5' and 3' regions から Kozak 配列など付加可能(任意)
4. Apply & Analyze をクリック

Review

Optimize (Optional)

Order

Example1

Add sequence ▾

Example1

Configure gene synthesis

Select vector and delivery quantity.

Note: Please check if insertion of kozak sequence is required for the vector chosen. You may refer to the kozak guide below the vector dropdown.

Missing a vector?

Vector: **pcDNA 3.1(+)**

View vector info | Cloning guide | Kozak guide

Insertion points: **HindIII [AAGCTT] - EcoRI [GAAT...]**

More info

Edit 5' and 3' regions

☐ Apply changes to all sequences ⓘ

Apply & Analyze

Sequence View

Plasmid View

Status ✓

Warning: You made changes to the configuration. When you're ready, clicking on 'Apply and Analyze' will apply them.

ATGGTGAGCAAGGGCGAGGAGCTGTTCACCGGGGTGGTCCCATCTCTGGTGGAGCTGGACGGCGACGTAACGGCCACAAGTTCAAGCGTGTCCGGCGAGGGCGAGGGCG
TACCACTCGTTCCTCGTCCGCTCCTCGACAAAGTGGGCCCCACCGGGTAGGACACGCTCGACCTGCCGTGCAATTTGCCGGTGTTCAGTGCACAGGCGCTCCCGCTCCCGC
ATGCCACCTACGGCAAGCTGACCTGAAGTTCATCTGCACACCGGCAAGCTGCCGTGCCCTGGCCCCACCTCTGTGACCACTTCACCTACGGCGTGCAGTCTTCGC
TACGGTGGATGCCGTTTCACTGGGACTTCAAGTAGACGTGGTGGCGCTTCGACGGGCACGGGACCGGGTGGGAGCACTGGTGAAGTGGATGCCGACGTACGGAAGCG
110 120 130 140 150 160 170 180 190 200 210
CCGCTACCCCGACCATGAAGCAGCAGCACTTCTTCAAGTCCGCCATGCCGAAGGCTACGTCCAGGAGCGCACCATCTTCTTCAAGGACGACGGCACTACAAGACC
GGCGATGGGGCTGGTGTACTTCGTCTGCTGAAGAAGTTCAGGCGGTACGGGCTTCGATGCAGGTCTCGCGTGGTGAAGAAGTTCCTGCTGCCGTTGATGTTCTGG
220 230 240 250 260 270 280 290 300 310 320
CGCGCCGAGGTGAAGTTCAGGGCGACACCTGGTGAACCGCATCGAGCTGAAGGGCATCGACTTCAAGGAGGACGGCAACATCTGGGGCACAAGCTGGAGTACAAC
GGCGGCTCCACTTCAAGCTCCCGCTGGGGACCACTTGGCGTAGCTCGACTTCCCGTAGCTGAAGTTCCTCTGCCGTTTGAAGACCCCGTGTTCGACCTCATGTTGA
330 340 350 360 370 380 390 400 410 420 430
ACAACGCCACAAGTCTATACCCGCCGACAAGCAAGAACGGCATCAAGTGAAGTTCAGAGCCGCCACAACATCGAGGACGGCAGCGTGCAGCTGCCGACCA
TGTGTGGTGTTCAGATATAGTGGCGCTGTCTGCTTCTTGGCGTAGTTCACCTTGAAGTTCAGGCGGTGTGTAGCTCCTGCCGTCGACGTGACGCGGTGGT
440 450 460 470 480 490 500 510 520 530 540
CTACCGCAGAACACCCCATCGGCGACGGCCCCGTGCTGCTGCCGACACCACTACCTGAGCACCCAGTCCGCCCTGAGCAAGACCCCAACGAGAAGCGCGATCAC
GATGGTCTGCTTGTGGGGGTAGCGCTGCCGGGGCAGCAGCGGGCTGTTGGTGATGGACTCGTGGGTGAGGGGAGCTGTTTCTGGGGTGTCTTTCGCGCTAGTG
550 560 570 580 590 600 610 620 630
BsrGI
ATGGTCTGCTGGAGTTCGTGACCGCCCGGGGATCACTCTCGGCATGGACGAGCTGTACAAGTAA
TACCAGGACGACCTCAAGCACTGGGGCGGCCCTAGTGAGAGCGCTACCTGCTCGACATGTTATT

Legend: X
EcoRI Cloning Site
Protected Region
EcoRI Unwanted Motif

6

研究用にのみ使用できます。診断用には使用いただけません。

Optimize (optional) 画面

1. 生物種を選択
2. ORF を選択
3. Optimize Sequence をクリック→ Success と表示されれば完了

Review

Optimize (optional)

Order

Example1

Add sequence ▾

Example1

Optimize sequence ⓘ

Optimize your sequence for better expression or to fix potential problems.

ⓘ Please select a Host Organism and ORF to optimize your sequence.

1

Host Organism
Homo sapiens ▾

2

ORF
1 - 720 ▾

Clear

Apply

Protect my cloning sites ⓘ

Protected Regions ⓘ

Motifs to Avoid ⓘ

3

Optimize Sequence

Sequence View

Plasmid View

Quality Graphs

Status ✓

✓ Success: Download and review cloned construct genbank file.

ATGGTGAGCAAGGGCGAGGAGCTGTTCACCGGGTGGTCCCATCTTGGTCGAGCTGGACGGCGACGTAACGGCCACAAGTTCAGCGTGTCCGGCGAGG
TACCACTCGTTCCCGCTCCTCGACAAGTGGCCCCACCAAGGAGGACGAGCTCGACCTGCCGCTGATTTGCCGTTTCAAGTCGACAGGCCGCTCC
M V S K G E E L F T G V V P I L V E L D G D V N G H K F S V S G E G
10 20 30 40 50 60 70 80 90 100
GCGAGGGCGATGCCACCTACGGCAAGCTGACCTGAAGTTCATCTGCACACCGGCAAGCTGCCGTCGCCCTGGCCACCCCTCGTGACCACTTCACCTA
CGCTCCCGCTACGGTGGATGCCGTTGACTGGGACTTCAAGTAGACGTGGTGGCGTTTCGACGGGCACGGGACCGGGTGGGAGCACTGGTGAAGTGGAT
G E G D A T Y G K L T L K F I C T T G K L P V P W P T L V T T F T Y
110 120 130 140 150 160 170 180 190 200
CGGCGTGAGTGTTCGCCCGCTACCCGACCATGAAGCAGCAGCACTTCTTCAAGTCCGCCATGCCCGAAGGCTACGTCCAGGAGCGCACCATCTTC
GCCGCACGTACGAAGCGGGGATGGGGCTGGTGTACTTCGTCTGTGTAAGAAGTTCAGGCGGTACGGGCTCCGATGCAGGTCTCGCGTGGTAGAAG
Y G V Q C F A R Y P D H M K Q H D F F K S A M P E G Y V Q E R T I F I
210 220 230 240 250 260 270 280 290 300
TTCAAGGACGACGCAACTACAAGACCGCGCCGAGGTGAAGTTCGAGGGCGACACCTGGTGAACCGCATCGAGCTGAAGGGCATCGACTTCAAGGAGG
AAGTTCCTGTGCGTTGATGTTCTGGGCGCGCTCCACTTCAAGCTCCCGCTGTGGGACCACTTGGCGTAGCTCGACTTCCCGTAGCTGAAGTTCCTCC
F K D D G N Y K T R A E V K F E G D T L V N R I E L K G I D F K E D
310 320 330 340 350 360 370 380 390 400
ACGGCAACATCCTGGGGACAAGCTGGAGTACAACACAGCCACAAGGTCTATATACCGCCGACAAGCAGAAGAAGCGCATCAAGGTGAACCTCAA
TGCCGTTGTAGGACCCCGTTCGACCTCATGTTGATGTTGCGGTGTTCCAGATATAGTGGCGGCTGTTGCTCTTTCGCGTAGTTCCTGAGTGAAGTT
D G N I L G H K L E Y N Y N S H K V Y I T A D K Q K N G I K V N F K

7 研究用にのみ使用できます。診断用には使用いただけません。

Order 画面

1. Choose add-on and fast delivery をクリック→Add-on service 画面が出る
2. Delivery Quantity にてプラスミド精製オプションを選択可能
(Add-on Service 画面にてその他のオプションの追加も可能)
3. Save をクリック

Review		Optimize (Optional)		Order
Your Sequence Configuration		Your Products ⓘ		
Add sequence ▾				
☐	Choose Add-on & Fast Delivery for Selected		Delete Selected	
☐	Example1 pcDNA 3.1(+) - Ampicillin 5µg dried plasmid DNA, Research Grade	720 bp	JPY 26,660 8 business days	① Choose add-on and fast delivery



Add-on Services

Choose the add-on services for your cloned gene.

Delivery Quantity

② TE, up to 100µg yield (midiprep) ▾

Production Time ⓘ

Standard ▾

☐ Sanger Sequencing ⓘ

☐ TSE-Free ⓘ

☐ Glycerol Stock ⓘ

☐ Stock Agar ⓘ

③

Cancel Save

Order 画面

1. Download Summary で合成配列やプラスミドマップの確認可能
2. Add to Cart から注文手続きへ

Review	Optimize (Optional)	Order				
Your Sequence Configuration Add sequence ▾		Your Products ⓘ				
☐ Choose Add-on & Fast Delivery for Selected Delete Selected						
☐ Example1 720 bp pcDNA 3.1(+) - Ampicillin TE, up to 100µg yield (midiprep), Transfection Grade		JPY 39,860 14 business days + Choose add-on and fast delivery				
		JPY 39,860.00 Project: Example1 Status: Draft <table><thead><tr><th>Item</th><th>Quantity</th></tr></thead><tbody><tr><td>Cloned genes</td><td>1</td></tr></tbody></table> Estimated production time ⓘ 14 to 14 business days Email will be sent when Quality Assurance files are available in Thermo Fisher Connect ⓘ Kindly review the plasmid sequence(s) before ordering ① Download Summary ▾ ② Add to Cart (1)	Item	Quantity	Cloned genes	1
Item	Quantity					
Cloned genes	1					

Thank you

研究用에만使用できます。診断용에는使用いただけません。

© 2024 Thermo Fisher Scientific Inc. All rights reserved.

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

実際の価格は、弊社販売代理店までお問い合わせください。

価格、製品の仕様、外観、記載内容は予告なしに変更する場合がありますのであらかじめご了承ください。

標準販売条件はこちらをご覧ください。 thermofisher.com/jp-tc

