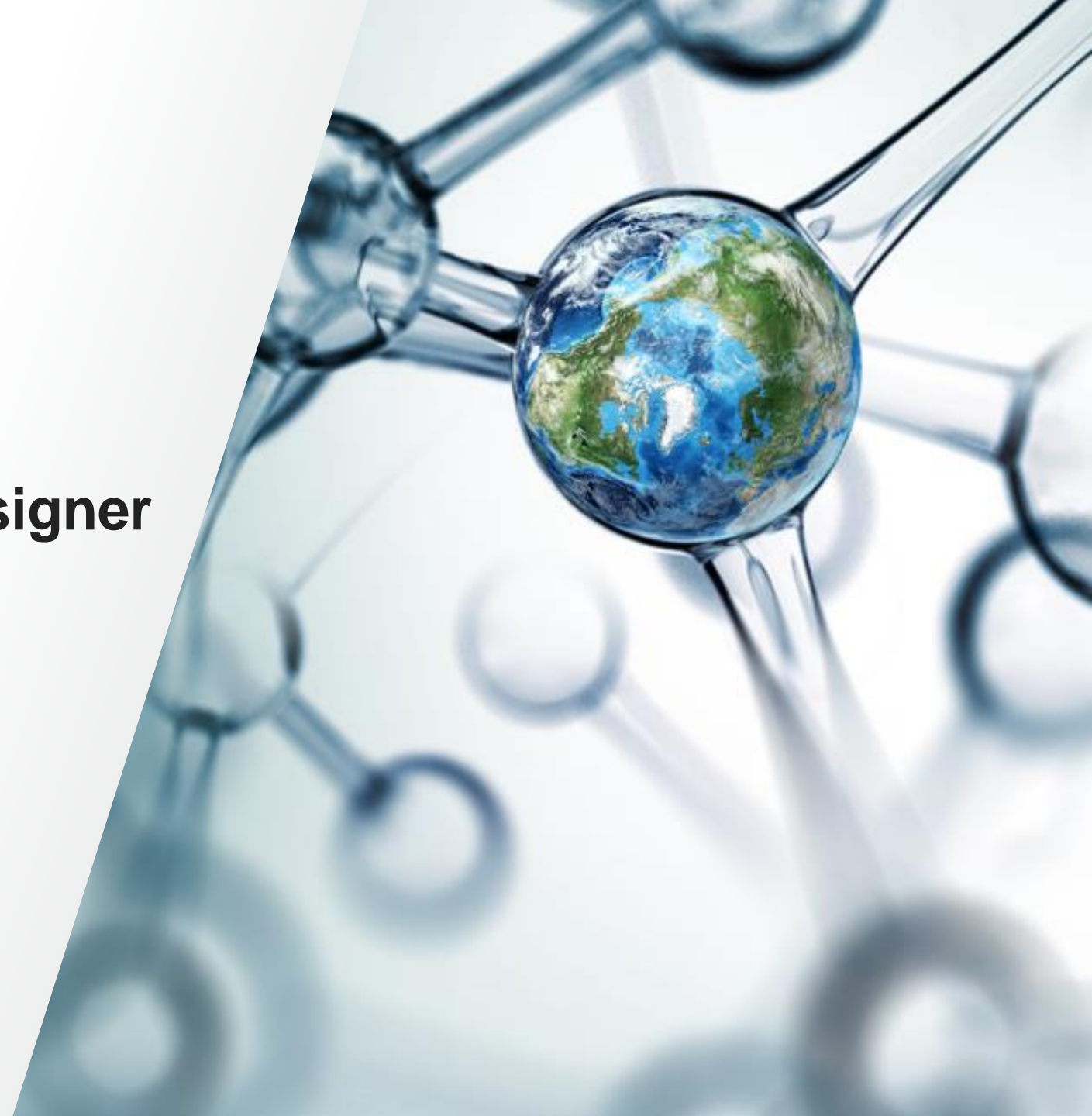


Invitrogen GeneArt Instant Designer オーダークイックガイド

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

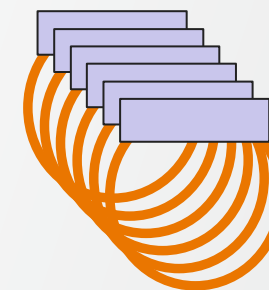
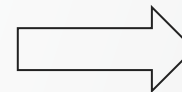
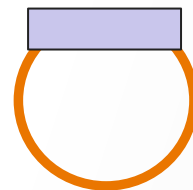
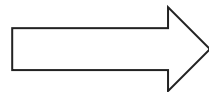
こちらは以下の内容の Web 入力クイックガイドです。(塩基配列の入力からプラスミド精製入力)

塩基配列入力

ベクター導入

プラスミド精製

ATGGTGAGC...



- 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。
- 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>
- 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

<input type="text" value="Example1"/>	<input type="text" value="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

<input type="checkbox"/> Sequence Name	DNA or Protein Sequence Important Information for Cloning and Protein Sequence(s)	5' Region (Optional)	3' Region (Optional)
<input type="checkbox"/> Example1	ATGGTGAGCAAGGGCGAGGAGCTGTTACCCGGGGTGGTCCCATCTGGTCGAGCTGGACGGCGACGTA...	Click to enter 5' region	Click to enter 3' region
<input type="checkbox"/>			
<input type="checkbox"/>			
<input type="checkbox"/>			
<input type="checkbox"/>			
<input type="checkbox"/>			

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.](#)

4

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. プラスミド精製オプションを追加可能
4. Edit 5' and 3' regions から Kozak 配列など付加可能(任意)
5. Apply & Analyze をクリック

The screenshot displays the software interface with three main tabs: Review, Optimize (optional), and Order. The Review tab is currently selected and shows the following configuration options:

- Example1**: Add sequence ▾
- Example1
- 1** Vector: **pcDNA 3.1(+)** (Missing a vector?)
- View vector info | Cloning guide
- 2** Insertion points: **HindIII [AAGCTT] - EcoRI [GAAT...]** (More info)
- 3** Delivery Quantity: **TE, up to 100µg yield (midiprep), ...**
- 4** Edit 5' and 3' regions
- Apply changes to all sequences ⓘ
- 5** **Apply & Analyze**

The Order tab is also visible, showing a **Warning**: "You made changes to the configuration. When you're ready, clicking on 'Apply and Analyze' will apply them." Below the warning is a sequence view with a plasmid map and sequence details.

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 i

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

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

Host Organism
Homo sapiens ▾

ORF
1 - 720 ▾

Clear
Apply

Protect my cloning sites i

Protected Regions i

Motifs to Avoid i

Optimize Sequence

Sequence View
Plasmid View
Quality Graphs
Status ✔

✔ Success: [Download](#) and review cloned construct genbank file.

```

ATGGTGAGCAAGGGCGAGGAGCTGTTCAACCGGGTGGTCCCATCTGGTTCGAGCTGGACGGCGACGTAACGGCCACAAGTTCAGCGTGTCCGGCGAGG
TACCACTCGTTCCTCCGCTCCTCGACAAGTGGCCCAACACGGGTAGGACAGCTCGACCTGCCGCTGCATTTGCCGGTGTCAAGTCGCACAGGCCGCTCC
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
GCGAGGGCGATGCCACCTACGGCAAGCTGACCCCTGAAGTTCATCTGCACACCGGCAAGCTGCCCGTCCCTGGCCACCCCTCGTGACCACCTTCACTTA
CGTCCCGCTACGGTGGATGCCGTTGACTGGGACTTCAAGTAGACGTGGTGGCCGTTTCGACGGGCACGGGACCGGGTGGAGCACTGGTGAAGTGGAT
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
CGGCGTGCAGTGTTCGCCCGCTACCCGACCATGAAGCAGCAGCACTTCTTCAAGTCCGCCATGCCCGAAGGCTACGTCAGGAGCGCACCATCTTC
GCCGACGTCACGAAGCGGGCGATGGGGCTGGTGTACTTCTGCTGTGCTGAAGAAGTTCAGGCGGTACGGGCTCCGATGCAGGTCCTCGCGTGGTAGAAG
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 F
210 220 230 240 250 260 270 280 290 300
TTCAAGGACGACGGCAACTACAAGACCCGCGCCGAGGTGAAGTTCGAGGGCGACACCTGGTGAACCGCATCGAGCTGAAGGGCATCGACTTCAAGGAGG
AAGTTCCTGCTGCCGTTGATGTTCTGGGCGCGCTCCACTTCAAGCTCCCGCTGTTGGGACCACTTGGCGTAGCTCGACTTCCCGTAGCTGAAGTTCCTCC
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
ACGGCAACATCCTGGGGCAAGCTGGAGTACAACACTACAACAGCCACAAGGTCTATATACCGCCGACAAGCAGAAGAACGGCATCAAGGTGAACCTCAA
TGCCGTTGATAGGACCCGTTTCGACCTCATGTTGATGTTGCGGTGTTCCAGATATAGTGGCGGCTGTTGCTCTTCCGCTAGTTCCTCAAGTTC
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

```

①

②

③

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		Your Products ⓘ	
Add sequence ▾			
<input type="checkbox"/> Choose Add-on & Fast Delivery for Selected Delete Selected			
<input type="checkbox"/>	Example1 pcDNA 3.1(+)-Ampicillin TE, up to 100µg yield (midiprep), Transfection Grade	720 bp	JPY 39,860 14 business days Choose add-on and fast delivery

JPY 39,860.00

Project: Example1
Status: Draft

Item	Quantity
Cloned genes	1

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

[Download Summary](#)

Add to Cart (1)

[Contact us](#)

①

②

Thank you

研究用にものみ使用できます。診断用には使用いただけません。

© 2023 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](https://www.thermofisher.com/jp-tc)

