

## Invitrogen<sup>™</sup> GeneArt<sup>™</sup> Instant Designerご利用マニュアル

バイオサイエンス事業部

2024年11月21日版

The world leader in serving science

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



**Thermo Fisher** S C I E N T I F I C

# **GeneArt DashBoard**

2 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

#### GeneArt Dashboard サインイン



- thermofisher.com/genesynthesis にアクセスし、[GeneArt Instant Designer で遺伝子合成をオーダー]の [注文する] をクリック
- 2 ユーザー名とパスワードを入力しサインイン
- 3 GeneArt Dashboardに移動
  - A) [Home]:各種サービスを選択(詳細は<u>P4</u>)
  - B) [Orders & Drafts]: プロジェクト概要を確認(詳細は<u>P41, P42, P43</u>)
  - c) [Project Templates]:複数配列一括設定ひな形を確認・修正(詳細はP23, P35)



2	3A			3B			
サインイン	GeneArt <sub>Welcome</sub>	New! Templates and bulk configuration features now make placing your larger o Watch for the option to configure and optimize multiple sequences during	frequent orders simple. X the import step.	GeneArt	Drafts 📾	Search by project name or proj	ect id Q
ユーザー名: *  次へ	Home     Grass & Grass     Fill     Project Templates     Support     Support     Priorig	Cloned Genes (100-12000 bp) (200-3000 bp)	Reorder Plasmid Order stored pasmid	Fiome     Orders & Drafts     Open Orders     Pact     Project Templates	Project Name         Project Id           1         Tiss1_Strings         2023AJUU0F           1         Tiss1_Strings         2023AJUU0F           1         D/CZapha         2023AJVED           1         2106 and Skip Sklus         2023AJVED           1         Manus Screenshot         2023AVED           1         Kau Skgenes         2023AVED           1         Gane Order contim         2023AVED	# Sequences         Project Type         Last modified         Project Type         Last modified         Project Type         Last modified         Project Type         Project Typ	Status 0 Drat A Drat Drat Drat Drat Drat Drat Drat Drat
サインインにお困りですか?		Start New Project Start New Project Learn more Learn more	Order Now Learn more	<ul> <li>Support</li> <li>Pricing</li> </ul>	Image: Constraint of Constraints         Constraints <thco< td=""><td>1         Cloned Genes         11-Aug-2023           3         Cloned Genes         03-Aug-2023</td><td>Dreft Dreft</td></thco<>	1         Cloned Genes         11-Aug-2023           3         Cloned Genes         03-Aug-2023	Dreft Dreft
		Custom Proteins Expressed and purified Directed E	Praries	GeneA 3 Celcome + Home	Add or edit project Templ Add or edit project templates w easy submission of future proje and select to configure multiple	ates - Cloned Genes th your favorite vector configurations and optimiza cts. To use a template, start a new cloned genes pr sequences during the import step.	3 tion settings to enable fast an oject from the Dashboard Ho
		Regest a Cable Learn more Learn	Guide	📜 Orders & Draf	fts  Actions Building Delete Configure	Template Name 🖕 Manual 230830	Last Mo 30-Aug-
関連ページ: <u>P4, P23, P35, P41, P42, P43</u>				H Project Tem Cloned Genes DNA Fragment	s		
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							

 $(\mathbf{1})$ 

### GeneArt Dashboard サービスの選択

- **Cloned Genes DNA Fragments Reorder Plasmid** (100-12000 bp) (200-3000 bp) Order stored plasmid Start New Project Start New Project Order Now Learn more Learn more Learn more AB $(\mathbf{C})$ Plasmid Reorder - Instant Designer 1. Enter project information (Optional) Default project name will be given if not specified Select the constructs for reordering Cancel Import & Save Project name Optional description onstructs that are not available for reorder are displayed in grey 😗 Constructs Oltems My Projects and Constructs Search by project name or id Search by construct name or id Q 2. Select sequence input method Download Excel template Construct Name Project Name Last Updated Construct Id No constructs selected Add → Upload Sequence File Manual Creatio ← Remove FASTA, GenBank or Excel format Copy & paste your sequences directly into Instant Designer
- GeneArt Dashboardでは次のサービスを選択可能
  - A) Cloned Genes→人工遺伝子合成(P6に続く)
  - B) DNA Fragments→GeneArt Strings DNA Fragments(P25Iに続く)
  - C) Reorder Plasmid→再注文、過去に合成した遺伝子のプラスミド精 製(P37に続く)
- 注意:Custom Proteins(タンパク質発現精製)、DNA Libraries(変異ライブラリー)はオンラインで対応しておりま せん。ご希望の場合はjpcustom@thermofisher.com(受 託サービス)までご連絡ください。担当者から連絡いたしま す。



#### 関連ページ:<u>P3, P6, P25, P37</u>

# GeneArt Instant Designer 人工遺伝子合成

## 新規プロジェクト設定画面 人工遺伝子合成

- プロジェクト名があれば [Project name] に、その他備考があれば[Optional description] に入力\*1(どちらも必須ではない)。前者が未入力の場合はProject ID (例:2023ABCDEF)が自動表示される
- 2 プロジェクトに関連タグを一つ付加することが可能。
- ③ [Download Excel template] をクリックして、配列を入力するファイルをダウンロード
- ④ ダウンロードしたファイルに以下の情報を入力。複数配列入力可能
  - A) Sequence Name(遺伝子名):半角英数字、記号は[-]と[\_]のみ、スペースは不可、20文字以内
  - B) DNA / Protein Sequence: DNA: ATGC表記、アミノ酸: 1文字表記のみ、アミノ酸配列の場合のストップコドンは アスタリスク【\*】
  - c) 5' Region (optional)(5'付加配列: Kozakなど): DNA: ATGC表記
  - D) 3' Region (optional)(3'付加配列:ストップコドンなど): DNA: ATGC表記

注意:クローニング用制限酵素サイトが入力・付加されていない場合は、後に選択した制限酵素サイトが 自動的に合成予定配列に付加される(P17参照)。配列中に存在する制限酵素サイトは利用できない。 最適化で合成予定配列中の制限酵素サイトの削除が可能(P10④)。

- 5 ファイルをアップロードするため、左下の [Upload Sequence File]をクリック
- 6 ファイルのアップロード画面に移動(<u>P7</u>に続く)
- ⑦ 配列情報を直接手入力する場合は右下の [Manual Creation] をクリック
- 8 ④と同様に配列情報を入力して右上の [Continue] をクリック(<u>P8</u>に続く)

\*1:全角文字・スペースはエラーの原因となりますので、ご注意ください。

関連ページ:<u>P4, P7, P8, P10, P17</u>

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





## 配列情報ファイルをアップロード 人工遺伝子合成

- 配列情報を入力したファイルをDrag、あるいは [Choose file] をクリックしてファイルを選択する と、自動的に配列情報がアップロードされる
- 2 アップロードが完了すると右側にファイルが表示される。ファイルは複数アップロード可能
- ③ アップロードが完了したら、右上の [Continue] をクリック
- ④ 入力情報が不十分な場合、[Check Status]を クリックして確認画面へと進む(<u>P8</u>に続く)
- 5 入力情報が充分な場合、[Stay on this page]を クリックして④と同じ確認画面へ進む(<u>P8</u>に続 く)、または[Continue to configure]をクリックし て[Review]へと進む。(<u>P9</u>に続く)



Check Status

**Thermo Fisher** 

SCIENTI

#### 関連ページ:<u>P6</u>, <u>P8</u>, <u>P9</u>

## 配列情報の確認・修正、最適化・ベクター・納期短縮の設定

- 入力した配列情報のリストが表示されるので、内容を確認。エラー(赤い[!])がある場合は[Configure]をクリック
- [Sequence Info] タブで配列情報の修正変更が可能
- ③ [Optimization Settings]で最適化の設定が可能。アミノ 酸配列は最適化が必須。最適化生物種のリストはP12, P13参照。
- (Clone configuration)タブ でベクターの種類・納期短縮オプションの選択が可能
- ⑤ エラーがないもの(緑の[✔])も変更可能。全て設定が 完了したら、右上の [Import & Save] をクリックして Review・Optimize・Order画面へ移動(P9へ続く)
- 6 複数遺伝子への一括設定適用は<u>P20</u>~参照
- プラスミド精製はP11, P18のOrderタブのAdd-on Servicesで設定可能参照(ここでは不可)

Project: 2023AAY2ND				
$\bigtriangledown$	2			
Upload Sequence(s)	Verify & Configure			<b>(5)</b>
2 Sequence(s) Errors (1)				Add Sequence - Import & Save
Configure Selected Delete Selected	Protein sequences must have a	host organism selected. You can do this fo	or one or multiple sequences.	
Actions Status	Sequence Name Vector 🛓	Delivery Quantity 🛓 Type	Host Organism 🛓	Length 🟮 DNA/Protein Sequence
🔲 🔟 Delete 🖉 Configure 🔗	HsActin_nuc Standard (pM	1X) 5μg dried plasmid DNA DNA	No optimization	1152 ATGGAATTCAAGCTTGG
Delete Configure	HsActin_aa Standard (pM	IX) 5μg dried plasmid DNA PRO.	No optimization	384 MEFKLGSCEEEDSTALV
Configure Sequence Sequence Info ▲ Optimization Settings Sequence name* Sequence type HsActin_aa F Region (Optiona) GAATTCGCCAAC GAATTCGCCAAC Sequence * MCEEEDSTALVCDNGSGLCKAGFAGDDAPRAVFPSINGRPRHK DEADSRAGILTLYYPIEHGIITNWDDMENWHHSTYNELFVAPEI ENMTGIMFETFNVPANYVAIDAVLSLVAGGRTTGIVLDSGDGVTT LDLAGRDLTDYLMRILTERGYSFVTTAEREIVRDIKEALCYVALDR Length:378	Clone configuration 、 ional) 記例情報ファイルを 2GVMVGMGGKDSYVG 当HPTLITEAPLINEKAINR HNPFYEGYALPHAIMR ENEMATAASSSSLEKS Cancel Save	Configure Sequence  Sequence Into  A Optimization Settings  Cone configurati Host Organism for Optimization  No optimization  A Press solid a North Optimization  Standard Motifs  Chick on the cloring gate to add or remove.  Standard Motifs  Chick o	Configure :     Sequence In     Sequence In     Sequence In     Over Please     chosen. You     Vector *     pcDNA3.4-TOPO     View vector info   Ck     Select insertion p     Production Time *     Standard     Add-on Servi     choose the add-on servi	Sequence ×  to Optimization Settings Clone configuration  c check if insertion of Kozak sequence is required for the vector × may refer to the kozak guide below the vector dropdown.  c consigning guide Kozak guide Missing a vector?  sints  c constraints  c cons
			CONTRACTOR OF CO	A · Sanger Sequencing () TSE-Free () Citycerol Stock () Stock Agar ()

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

8

関連ページ: P6, P7, P9, P11, P12, P13, P15, P18, P20, P21

#### Thermo Fisher

## Review ベクター・付加配列の編集

- 事前に選択したベクターが表示される。変更する場合は [Vector]から再度選択(詳細は<u>P14</u>, <u>P15</u>, <u>P16</u>参照)。制限酵 素法の場合は挿入部位を選択(詳細は<u>P17</u>参照)。
- [Edit 5' and 3' regions] をクリックすると5'付加配列、3'付加 配列の入力が可能(Kozak、制限酵素サイト、クローニング用 付加配列など)

注意:クローニング用制限酵素サイトが入力・付加されていない場合は、後 に選択した制限酵素サイトが自動的に合成予定配列に付加される(P17参 照)。配列中に存在する制限酵素サイトは利用できない。最適化で合成予 定配列中の制限酵素サイトの削除が可能(P10④)。

- ③ [Apply changes to all sequences] に☑を入れると、同一プロジェクト内の全ての遺伝子に、チェック画面のベクター・5'/3'付加配列の設定が適用される。付加配列が空欄の場合は全て空欄になるので注意。
- ④ [Apply and Analyze] をクリックして変更を反映
- ⑤ プラスミド精製はP11, P18のOrderタブのAdd-on Servicesで設定可能参照(ここでは不可)
- 6 塩基配列を直接修正する場合は該当部位を選択後に右上の [Edit]をクリック、枠内を修正後 [Apply] をクリック
- アミノ酸配列が表示されている場合は、該当アミノ酸をクリック。[Edit]をクリックすると塩基の置換が可能、枠内を変更後 [Apply]をクリック
- 8 [Plasmid View] でマップの確認が可能
- 9 複数遺伝子への一括設定適用は<u>P20</u>~参照

関連ページ: <u>P7, P8, P10, P11, P14, P15, P16, P17, P18, P20, P21</u>、<u>P22</u>

Review Optimize (Optional) Order (8) Plasmid View Status 🌄 Sequence View Test01 Configure gene synthesis Add sequence Select vector and delivery quantity. Warning: You made changes to the co icking on 'Apply and Analyze' will apply them Note: Please check if HsActin nuc insertion of kozak sequence is EcoRI required for the vector GAATTCGCCACCATGTGTGAAGAAGAGGACAGCAC TGGTGTGTGACAATGGCTCTGGGCTCTGTAAGGCCGG chosen. You may refer to the HsActin aa kozak guide below the vector dropdown. 10 40 RT TCCCATCCATTGTGGGACGTCCCAGACATCAGGGGGTGATGC ATGATGATGATGTGTG Missing a vector' 6 AGGGTAGGTAACACCCTGCAGGGTCTGTAGTCCCCCACTACC TCGCCACCATGCACCACCAC pcDNA3.4-TOPO Edit ACTACTACTACACAC CTTAAGCGGTGGTACGTGGTGGTGGT 150 View vector info Cloning guide Kozak guid 5'region GAAGTACCCGATAGAACATGGCATCATCACCAACTGGGACGAC CTTCATGGGCTATCTTGTACCGTAGTAGTGGTTGACCCTGCT 228 230 749 250 268 CACCACCACCACCACCAC Edit 5' and 3' regions CTGCTCACGGAGGCACCCCTGAACCCCAAGGCCAACCGGGAGA GACGAGTGCCTCCGTGGGGACTTGGGGTTCCGGTTGGCCCTC 5' Region (Optional) Length: 18 Start: 16 End: 33 Position: 34 GAATTCGCCACC 330 350 TGTCTCTCTATGCCTCTGGACGCACAACTGGCATCGTGCTGG/ 3' Region (Optional) ACAGAGAGATACGGAGACCTGCGTGTTGACCGTAGCACGACC 0 Cancel Apply AAGCTT 458 478 TTALLUTGTTTTTTTTTTTGTUGATGLALLUAUTGUT BglII GCGTCTGGATCTGGCTGGCCGAGATCTCACTGACTACCTCATG Apply changes to all sequences CGCAGACCTAGACCGACCGGCTCTAGAGTGACTGATGGAGTA Apply & Analyze 550 570 GCCACCATC LI I AAGCGGTGGTACGTGGTGGTGGTAGTGGTAACGCTCCTTCTCCT Length: 1152 Start: 1 End: 1 Selected: 0 5'region ⑤プラスミド精製はOrderタブのAdd-on Servicesで設定可能 CAT CATCG Add-on Services X GTAGC End: 18 Position: 19 Choose the add-on services for your cloned gene Start: 16 Length: 3 )I) Delivery Ouantit 5µg dried plasmid DN/ Sanger Sequencing ด Cancel Apply TSE-Free Glycerol Stock Standard Stock Agar 6

Cancel

Save

9 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 合成予定配列の最適化

- ① [Optimize] をクリック
- [Host Organism] に事前に選択したコドン最適化用生物種が 表示(変更可能、最適化生物種のリストはP12, P13参照)
- [ORF] をクリック、[Define manually] で最適化を行う範囲を 入力(3塩基単位)、または選択肢から指定
- ④ 3種類の条件を追加可能
  - Protect my cloning sites]:最適化時に変更したくない制限酵素サイトを指定可能
  - Protected Regions]:最適化時に変更したくない配列 領域を指定可能
  - [Motifs to Avoid]:最適化後の配列に含めたくない制
     限酵素サイトを指定可能、<u>遺伝子内の制限酵素サイト</u>
     をクローニングに利用する場合、この指定は必須(A)
- 5 [Optimize Sequence] をクリックして最適化を実行
- 6 [Plasmid View] でマップを確認可能
- ⑦ [Quality Graphs] で最適化後の結果を確認可能
- 8 複数遺伝子への一括設定適用は<u>P20</u>~参照



関連ページ:<u>P6, P9, P12, P13, P20, P21</u>

### Order プラスミド精製・オプション追加 カートに追加

- (1) [Order] をクリック
- 2 価格と作業期間を確認
- ③ チェックを入れて[Choose Add-on,,,]をクリックすると納 期短縮サービス(Express: 1日短縮、Super Speed: 2~ 5日短縮)、プラスミド精製(P18参照)、およびサンガー シーケンス・TSE-Free・大腸菌ストックのオプションが追 加可能。
- ④ 個々の遺伝子に対してオプションを追加する場合は [Choose Add-on,,,]をクリック
- ⑤ 必要に応じて [Download Summary] から配列情報およびレポートをダウンロード
- ⑥ [Add to Cart] をクリックして注文手続きへと進む(詳細は [GeneArt人工遺伝子合成サービス\_ご注文マニュアル] のP8[6 カート内での価格・予想処理時間の確認]参照)
- ⑦ 複数遺伝子への一括設定適用は<u>P20</u>~参照

Review Optimize (options		er				
Your Sequence Configuration	Your Products 0		JPY 139,264.00			
Choose Add-on & Fast Delivery for Selected Delete Selected	2		Status: Draft			
♥	JPY 62,256 14 business days Choose add-on and fast delive	ED/	Item Cloned genes Estimated production time () 8 to 14 business days	Quantity 3		
pcDNA 3.1(+) - Ampicillin 5µg dried plasmid DNA, Research Grade	JPY 38,420 8 business days Choose add-on and fast delive	ny	Email will be sent when Quality Assurar are available in Thermo Fisher Connect	uence(s)		
Image: marked state state     1146 bp       pcDNA 3.1(+) - Ampicillin     5µg dried plasmid DNA, Research Grade	JPY 38,588 8 business days Choose add-on and fast delive	ny	5 before ordering Download Summary 6 Add to Cart (5)		Download Summa Download excel fo Download report +	r <b>y</b> rmat - FASTA format
3	(	4				
Bulk Add-on Services Important: Fields that cannot be edited in bulk are not displayed. Changes will be applied to the selected sequences.	×	Add-on Services Choose the add-on services for	for your cloned gene.		×	
Choose the add-on services for your cloned gene. Delivery Quantity Syg dried plasmid DNA		Delivery Quantity 5µg dried plasmid DNA Production Time 0 Standard	Sanger Sequen     TSE-Free     Glycerol Stock     Stock Agar	cing () ()		
Cancel	Save			Cancel	Save	

#### 関連ページ:<u>P18</u>, <u>P20</u>, <u>P21</u>

#### 最適化生物種

 GeneArt Instant Designerで選択できる生物種は 右の通り

- ② リストにある生物種の学名と一般名は次ページを 参照
- ③ リストにない生物種での最適化は
   jpgene@thermofisher.com
   までお問い合わせください

search No optimization Arabidopsis thaliana Aspergillus niger Aspergillus oryzae Bacillus subtilis Bos taurus Brassica napus Caenorhabditis elegans Canis familiaris Clostridium acetobutylicum Corynebacterium glutamicum ATCC 13032 Cricetulus griseus Danio rerio Drosophila melanogaster Escherichia coli Gallus gallus

No optimization

#### Glycine max

.

Q

Homo sapiens

Hordeum vulgare

Hypocrea jecorina (Trichoderma reesei)

Kluyveromyces lactis 0406

Lactobacillus casei ATCC 334

Lactococcus lactis

Lycopersicon esculentum

Macaca mulatta

Mus musculus

Nicotiana benthamiana

Nicotiana tabacum

Oryctolagus cuniculus

Oryza sativa

Pichia angusta (Hansenula polymorpha)

Pichia pastoris

Plasmodium falciparum

Rattus norvegicus

Saccharomyces cerevisiae

Schizosaccharomyces pombe

Spodoptera frugiperda

Sus scrofa

Synechococcus elongatus

Trichoplusia ni

Triticum aestivum

Vaccinia virus

Xenopus laevis

Yarrowia lipolytica

Zea mays

関連ページ: <u>P8</u>, <u>P10</u>, <u>P20</u>, <u>P21</u>, <u>P27</u>, <u>P29</u>, <u>P32</u>, <u>P33</u>

#### 最適化生物種リスト

学名	一般名
No optimization	最適化なし
Arabidopsis thaliana	シロイヌナズナ
Aspergillus niger	黒色麹菌、黒麹カビ、クロコウジカビ
Aspergillus oryzae	日本麹カビ、ニホンコウジカビ
Bacillus subtilis	枯草菌
Bos taurus	牛、ウシ
Brassica napus	セイヨウアブラナ
Caenorhabditis elegans	線虫(C.elegans)
Canis familiaris	犬、イヌ
Clostridium acetobutylicum	クロストリジウム・アセトブチリクム
Corynebacterium glutamicum ATCC 13032	コリネバクテリウム・グルタミクム ATCC13032株
Cricetulus griseus	チャイニーズハムスター、CHO細胞
Danio rerio	ゼブラフィッシュ
Drosophila melanogaster	キイロショウジョウバエ
Escherichia coli	大腸菌
Gallus gallus	鶏、ニワトリ
Glycine max	大豆
Homo sapiens	人、ヒト
Hordeum vulgare	大麦
Hypocrea jecorina (Trichoderma reesei)	トリコデルマ・リーゼイ
Kluyveromyces lactis 0406	クルイウェロマイセス酵母
Lactobacillus casei ATCC 334	ラクトバチルス・カゼイ乳酸菌 ATCC 334株
Lactococcus lactis	ラクトコッカス・ラクティス、ラクチス乳酸菌、ラクチス菌

学名	一般名
Lycopersicon esculentum	トイト
Macaca mulatta	赤毛猿、アカゲザル
Mus musculs	ネズミ、マウス
Nicotiana benthamiana	ベンサミアナタバコ
Nicotiana tabacum	タバコ
Olyctolagus cuniculus	ウサギ
Oryza sativa	稲、イネ
Pichia angusta (Hansenula	ピキマ群母の一番
polymorpha)	して、時時の一種
Pichia pastoris	ピキア酵母の一種
Plasmodium falciparum	熱帯マラリア原虫
Rattus norvegicus	ラット
Saccharomyces cerevisiae	出芽酵母
Schizosaccharomyces pombe	分裂酵母
Spodoptera frugiperda	3トウガ、Sf9細胞、Sf21細胞
Sus scrofa	イノシシ・豚・ブタ
Synechococcus elongatus	シアノバクテリア(Engineering Kits, A14259)
Trichoplusia ni	イラクサギンウワバ、High Five™細胞
Triticum aestivum	パン小麦
Vaccinia virus	ワクシニアウイルス
Xenopus laevis	アフリカツメガエル
Yarrowia lipolytica	ヤロウイア酵母
Zea mays	トウモロコシ

関連ページ: <u>P8</u>, <u>P10</u>, <u>P20</u>, <u>P21</u>, <u>P27</u>, <u>P29</u>, <u>P32</u>, <u>P33</u>

#### 選択可能なベクターのリスト

 GeneArt Instant Designerで選択可能な ベクターは右の通り。以下の4種類に分類

- A) 標準ベクター(赤枠:pMX,pMA,pMK)
- B) Express cloningベクター(青枠:9種類の Invitrogenベクター)
- c) 13種類のInvitrogenベクター(緑枠)
- D) お客さまご提供のカスタムベクター(オレンジ 枠、追加方法はP15, P16参照)

※pMXベクターは、pUCベクターをベースにしたク ローニングベクターで発現ベクターではありません。 約95%のオーダーはアンピシリンまたはカナマイシ ン耐性遺伝子が組み込まれたベクターで納品されま すが、その他の薬剤耐性遺伝子が用いられる可能 性もあります。pMAやpMKを選択することにより、 マーカーを固定することも可能です(追加料金が必 要となります)。

※ Express cloningベクターへのExpress cloning は5 Kb以下の遺伝子が対象となります。なお pcDNA3.1(+)など制限酵素クローニングベクターは 5 Kb以上の遺伝子もサブクローニングでご注文可 能です。

関連ページ:<u>P8, P9, P15</u>, <u>P16, P20, P21</u>



#### 新規サブクローニングベクターの追加(1)

 $(\mathbf{3})\mathbf{A}$ 

- 1 新規ご提供ベクターをリストに追加する場合は、ベク ターリストの最下部にある[+ I want to add new vector(s)]をクリック
- 2 ベクター配列情報をアップデートする方法を次の2種類 から選択
- 3 [Upload Vector Sequence...]: ファイル(FASTA/ GenBank / Excel)をドラッグアンドドロップしてアップ ロード。アップロード用エクセルフォームをダウンロード 可能(A)。
- (4) [Copy & Paste Vector...]:ベクター名・全長塩基配列を 直接入力\*1
- (5) [Continue]をクリック、P16へ移動

\*1:全角文字・スペースはエラーの原因となりますので、ご注意ください。





#### 関連ページ: P9, P14, P16

## 新規サブクローニングベクターの追加(2)

- ① アップロードしたベクター配列情報をリストで表示
- [Edit]でベクター名を変更可能
- ③ [Plasmid View]でマップを確認可能
- ④ [Status]が緑の✔ならOK、ErrorやWarningの場合は次の点を満たしているかどうか確認
  - A) ベクター配列は薬剤耐性マーカーを含む
  - B) ベクター配列は複製起点Oriを含む
  - c) ベクター配列はATGCの4種のみ
  - D) ベクター名は半角英数字、記号は[-]と[\_]のみ
  - E) ベクター名にスペース・特殊な文字は含まれていない
  - F) ベクター名は15文字以内
- ⑤ [Continue to Cloned Genes]をクリック(P9へ続く)
- 6 ベクターリスト下部に追加
- ご注文後、弊社からMaterial Transfer Formを送付
- 8 ベクターとご入力・ご署名済みMaterial Transfer Form を弊社羽田事業所へご送付(詳細は<u>マニュアル[Option</u> <u>3 合成・構築・登録済みプラスミドのご利用方法、およ</u> びプラスミドのご提供方法]P13~16参照)



関連ページ:<u>P9</u>, <u>P14</u>, <u>P15</u>

#### Thermo Fisher

## サブクローニング用制限酵素サイトの選択

- [Insertion points] をクリッ クして、[Customize manually]を選択
- ② [Cloning Configuration]
   でクローニング用制限酵
   素サイトを選択
- ③ 表示されたマップで正しく クローニングされているこ とを確認(A)。制限酵素サ イトの再変更可能(B)
- ベクターとインサートの組 み合わせが正しくない場 合は右下の [other possible cloning strategies] をクリックする と他の候補が表示
- ⑤ 正しいマップを選択し、
   [Save] をクリック(P9に続く)



PciI [ACATGT]

Х



関連ページ:<u>P6</u>, <u>P9</u>, <u>P20</u>, <u>P21</u>,

### 選択可能なプラスミド精製 Add-on Servicesで設定可能

- 1 最少DNAは凍結乾燥DNA 5 µg
- ② プラスミド精製のスケールは以下の5種類から選択可能。 収量は目安で保証はなし。取得分を納品。培養スケー ルは固定
  - A) Midiprep (~100  $\mu$ g, 50 mL $\pi$  $\tau$ - $\mu$ )
  - B) Maxiprep (~500 μg, 250 mLスケール)
  - C) Megaprep (~1 mg, 500 mL $\lambda \tau$ - $\mu$ )

  - E) 2xGigaprep(10~15 mg, 5 Lスケール)
     精製スケール(収量の目安, 培養スケール)
- バッファーはTE, H2O, PBSの3種類から選択可能
- 4 Add-on Servicesで設定可能

Choose the add-on services for your doned gen Delivery Quantity  Sug dried plasmid DNA  Production Time  Standard	e. Sanger Sequencing () 1 TSE-Free () Glycerol Stock ()
---	--

能。	Delivery Quantity *	
	5µg lyophilized plasmid DNA, Research Grade	•
	search	Q
2	TE, up to 100µg yield (midiprep), Transfection Grade	•
3	TE, up to 500µg yield (maxiprep), Transfection Grade	
TE	TE, 1mg yield (megaprep), Transfection Grade	
	TE, 1-10 mg yield (gigaprep), Transfection Grade	
	TE, 10-15 mg yield (2x gigaprep), Transfection Grade	
2	H2O, up to 100µg yield (midiprep), Transfection Grade	PBS, up to 100µg yield (midiprep), Transfection Grade
3	H2O, up to 500µg yield (maxiprep), Transfection Grade	PBS, up to 500µg yield (maxiprep), Transfection Grade
H2O	H2O, 1mg yield (megaprep), Transfection Grade	PBS, 1mg yield (megaprep), Transfection Grade
	H2O, 1-10 mg yield (gigaprep), Transfection Grade	PBS, 1-10 mg yield (gigaprep), Transfection Grade
	H2O, 10-15 mg yield (2x gigaprep), Transfection Grade	PBS, 10-15 mg yield (2x gigaprep), Transfection Grade

#### 関連ページ: <u>P8</u>, <u>P9</u>, <u>P11</u>, <u>P20</u>, <u>P21</u>, <u>P38</u>

# 複数遺伝子への一括設定適用・設定の保存 人工遺伝子合成編

## 複数遺伝子への一括設定適用

#### **Thermo Fisher** S C | E N T | F | C

- 1番上のチェックボックスに☑を入れて、複数遺伝子を 一括選択
- ② [Configure Selected] をクリック
- ③ 二通りの選択肢から選択
  - A) [Configure sequences manually]:新規に設定
  - B) [Apply project template]: 既存の雛形から選択
- ④ [Continue]をクリック
- 5 設定画面へ移行。次の3つのタブで各種設定が可能
- 6 [Cloned Gene Settings]:
  - A) ベクター
  - B) クローニング用制限酵素サイト

C) 5'/3'付加配列

- ⑦ [Optimization Settings]
  - A) 最適化生物種
  - B) 除外する配列
- 8 [Add-ons]
  - A) プラスミド精製
  - B) 納期短縮オプション
  - c) 各種オプション(サンガー法・大腸菌グリセロールストック)

Manual 230830

関連ページ: P8, P9, P10, P11, P12, P13, P14, P17, P18



#### **Thermo Fisher** SCIENTI

### 各種設定・雛形として保存

	Cloned Gene Settings	Optimization Setti	ngs Add-ons				
	Vector 1	Missing a vector?	Insertion Points				
	Select vector	-	Select insertion point -				
	Cloning guide						
3	5' Region (Optional) 🚯		3' Region (Optional) 🚯				
	Enter 5' region		Enter 3' region				
		(	or)				
	Choose to delete any existing 5' & 3' region       Delete 5' Region       Delete 3' Region						
	Delete 5' Region	Delete 3' Regi	on				

- [Vector]: クローニング用ベクター (1)
- [Insertion Points]: クローニング用制限酵素サイ 2 F

3 [5'/3' Region (Optional)]:付加配列

	Cloned Gene Settings Optimization Setting	ngs Add-ons		Cloned Gene Settings
1	Host Organism for Optimization 0	1		Choose the add-on serv
	Homo sapiens -			Delivery Quantity
(2)	Dptimize the sequence upon Import & Save	0		TE, 1mg yield (megapi
				Production Time 🔒
(3)	Motifs to Avoid		2	Standard
	Define motifs to avoid in your sequences. Selected	I sites will be excluded during optimization.		Sanger Sequencing
	Standard Motifs Custom Motifs			4 TSE-Free 0
				5 Glycerol Stock 0
	Click on the cloning site to add or remove.			6 Stock Agar 🕕
	Standard Motifs	Standard Motifs To Avoid		<u> </u>
	Hind Q	Q	$\overline{7}$	Save settings as project
		Remult (COATCO)		Note: Vector, host organism,
	Tillian (AAGOTT)	BanningGokrooj		template for future projects.
		EcoRI [GAALIC]		Project Template Name *
			8	Manual 241114
(1)	[Host Organism]: 最近	<b>箇化生物</b> 種を選択。最		
	適化しない場合は[No o	ptimization]を選択		
(2)	Optimize the sequence	e 1:1を入れると	l	
	Import & Save 由に是近	高化を宇体		1 [Delivery (
		ᄢᇆᆁᅕᄴ ᇦᇆᆁᅕᄴ		(Production)
3	[IVIOUITS to Avoid]: 最適化	L 配列から 除外する 配列		3 [Sanger S
				(TSE-Free
				5 [Glycerol



- [Project Template Name]: 雛形の名前を入力\*1 (8)
- **(9**) [Save changes]: クリックして一括適用・雛形保存
- \*1:全角文字・スペースはエラーの原因となりますので、ご注意ください。

関連ページ: P8, P9, P10, P11, P12, P13, P14, P17, P18

### 複数遺伝子への一括設定適用の完了

- [Import & Save]をクリック、Reviewへ移動、この間に最 適化を実行(<u>P9</u>へ続く)

ſ	0.5								Add Sequence	Import & Save
	2 Seque	Figure Colorted III Dolote		Durbin erennen			- 46 - 6			
Ļ	✓ Con	ingure Selected 🔟 Delete		Protein sequence	s must have a host o	organism selected. You can d	o this for one	e or multiple sequences. 🕡		
•	~	Actions	Status	Sequence Name	Vector 👙	Delivery Quantity 🍦	Туре 🌲	Host Organism 🍦	Length 🖯	DNA/Protein Sequence
	<b>_</b>	🛍 Delete 🖉 Configure	0	HsActin_nuc	pcDNA 3.1(+)	TE, 1mg yield (megapr	DNA	Homo sapiens	1152	ATGGAATTCAAGCTTGG
	<b>~</b>	🛍 Delete 🖉 Configure	0	HsActin_aa	pcDNA 3.1(+)	TE, 1mg yield (megapr	PRO	Homo sapiens	384	MEFKLGSCEEEDSTALV

#### 関連ページ:<u>P9</u>

### ー括設定雛形の確認・修正

Thermo Fisher

- ① [Project Templates]から[Cloned Genes] を選択
- ② [Template Name]から該当の雛形をクリックして内容を 確認
- ③ 修正する場合は[Configure]または[Edit]をクリック
- ④ 新規に作成する場合は[Add New Template]をクリック

関連ページ	:	<u>P3</u>
-------	---	-----------

GeneArt	Project Templat	tes - Cloned Ge	nes O	Add New Template
Welcome	Add or edit project templates with yo easy submission of future projects. T and select to configure multiple sequ	our favorite vector configurations and io use a template, start a new cloned uences during the import step.	l optimization settings to enable fas I genes project from the Dashboard	st and I Home
(1)	Actions 3	Template Name 🌲	Las	t Modified 🖕
Orders & Drafts	Delete 🖉 Configure	Manual 230830	30-/	Aug-2023
Cloned Genes DNA Fragments				
GeneArt Welcome	<sup>₄ Back</sup> Manual24	41114	3	) Edit
A Home	Cloned Gene Set	tings Optimization Settin	ngs Add-ons	
🗮 Orders & Drafts	Vector:		Insertion Points:	
# Project Templates	5' Region:		3' Region:	
Cloned Genes DNA Fragments	-		- /	
GeneArt Edit	Project Template	emplate that can be applied to future Cloned Genes	projects.	
Home     Manual	mplate Name* 241114			
Torders & Drafts Cloned	d Gene Settings Optimization Settings Add-	ons		
H Project Templates  Cloned Genes	Note: Please check if insertion of kozak sequence is re	equired for the vector chosen. You may refer to the k	sozak guide below the vector dropdown. $ imes$	
File Exchange Vector	Mise	sing a vector? Insertion Points		
PcDI	NA3.4-TOPO	Select insertion point	*	
Pricing 5' Denii	guide   Kozak guide	3' Beninn (Ontional) 🚔		
- 5 Regit	r 5' region	Enter 3' region		
	-			

# **GeneArt Strings DNA Fragments**

#### **Thermo Fisher** S C I E N T I F I C

### GeneArt Strings DNA Fragments 新規プロジェクト設定画面

- プロジェクト名があれば [Project name] に、その他備考があれば [Optional Description] に入力\*1(どちらも必須ではない)。前者が 未入力の場合はProject ID(例:2023ABCDEF)が自動表示される
- 2 プロジェクトに関連タグを一つ付加することが可能。
- ③ [Download Excel template] をクリックして、配列を入力するファイ ルをダウンロード
- ④ ダウンロードしたファイルに以下の情報を入力。複数配列入力可能
  - A) Sequence Name(遺伝子名):半角英数字、記号は[-]と[\_]のみ、スペース は不可、20文字以内
  - B) DNA / Protein Sequence: DNA: ATGC表記、アミノ酸: 1文字表記のみ、アミノ酸配列の場合のストップコドンはアスタリスク【\*】
- 5 ファイルをアップロードするため、左下の [Upload Sequence File] をクリック
- 6 ファイルのアップロード画面に移動(P26に続く)
- ⑦ 配列情報を直接手入力する場合は右下の [Copy and paste] をク リック
- ⑧ ④と同様に配列情報を入力して右上の [Continue] をクリック(P27)
   に続く)

\*1:全角文字・スペースはエラーの原因となりますので、ご注意ください。

関連ページ:<u>P4</u>, <u>P26</u>, <u>P27</u>





#### GeneArt Strings DNA Fragments 配列情報ファイルをアップロード

- 配列情報を入力したファイルをDrag、あるいは [Choose file] をクリックしてファイルを選択する と、自動的に配列情報がアップロードされる
- ② アップロードが完了すると右側にファイルが表示される。ファイルは複数アップロード可能
- ③ アップロードが完了したら、右上の [Continue] をクリック
- ④ 入力情報が不十分な場合、[Check Status]を クリックして確認画面へと進む(<u>P27</u>に続く)
- 5 入力情報が充分な場合、[Stay on this page]を クリックして④と同じ確認画面へ進む(<u>P27</u>に続 く)、または[Continue to configure]をクリックし て[Review]へと進む。(<u>P28</u>に続く)



**Thermo Fisher** 

SCIENTI

関連ページ:<u>P25</u>, <u>P27</u>, <u>P28</u>

### GeneArt Strings DNA Fragments 配列情報の確認・修正

- 入力した配列情報のリストが表示されるので、内容を確認。
   エラー(赤い[!])がある場合は[Configure]をクリック(A)
- [Sequence Info] タブで配列情報の変更が可能。③のよう にエラーメッセージのある項目は修正が必須。
- 3 アミノ酸配列は最適化が必須。最適化生物種のリストは P12, P13参照。
- ④ エラーがないもの(緑の[✔])も修正変更可能。全て設定が 完了したら、右上の [Import & Save] をクリックしてOrder 画面へ移動(P28)
- 5 複数配列を一括設定する場合はP32~を参照

	Project: Strings_Manual_230831		
Ŭ	Upload Sequences	2 Verify & Configure	
,	4 Sequence(s) Errors (2)		
0	🖉 Configure Selected 🔟 Delete Selected	Protein sequence(s) must have a hos	t organism selected. You can do this for one or <b>multiple sequences ()</b>
	Actions (A) Status	Sequence Name Type +	ength 1 DNA/Protein Sequence
	📄 🗑 Delete 🖉 Configure 🥥	HsActin_nuc DNA 1	152 ATGGAATTCAAGCTTGGATCCTGTGAAGAAGAGGACAGCACTGCCT
,	🗌 🗑 Delete 🖉 Configure 🏮	HsActin_aa PROTEIN 3	84 MEFKLGSCEEEDSTALVCDNGSGLCKAGFAGDDAPRAVFPSIVGRPR
	🔲 🖻 Delete 🖉 Configure 🥥	HsActin_nuc DNA 1	152 ATGGAATTCAAGCTTGGATCCTGTGAAGAAGAGGACAGCACTGCCT
	📄 🗑 Delete 🖉 Configure 🌒	HsActin_aa PROTEIN 34	84 MEFKLGSCEEEDSTALVCDNGSGLCKAGFAGDDAPRAVFPSIVGRPR
ر 3 (3	Sequence info Sequence name * HsActin_aa 5' Region(Optional) GCCACC Host Organism * No optimization Required field Sequence * MEFKLGSCEEEDSTALVCDNC VGMGQKDSYVGDEAQSKRG PEEHPTLITEAPLNPKANREKL LDSGDGVTHNVPIYEGYALPH EIVRDIKEKLCYVALDFENEMA Length: 384	Sequence type * PROTEIN 3' Region(Optional) TAG Host Organism * No optimization Host Organisms Homo sapiens Hordeum vulgare	



Import & Save

**Thermo Fisher** 

SCIENT

Add Sequence +

(**4**)

Import & Save

Host Organism

No optimization

No optimization

Add Sequence -

### GeneArt Strings DNA Fragments Order 注文可否の確認

- 入力した配列情報がGeneArt Strings DNA Fragments で注文可能か確認。価格・作業期間が表示されている ものは可能
- ② [Sequence cannot,,,]は注文不可。次の二通りの選択 が可能
  - A) [Fix/optimize to get string]:最適化・配列修正を実施(P29へ続く)
  - B) [Order as cloned gene]:人工遺伝子合成(pMXベクター・5 µgに限定)に変更(P30へ続く)



**Thermo Fisher** 

SCIENTI

関連ページ: <u>P26</u>, <u>P29</u>, <u>P30</u>

### GeneArt Strings DNA Fragments Optimize 最適化

- [Show problems] をクリックして、問題部位を表示。右配列中の赤矢印が問題部位(A)
- [Host Organism] に事前に選択したコドン最適化用生物種が 表示(変更可能、最適化生物種のリストは<u>P12</u>, P13参照)
- ③ [ORF] をクリック、[Define ORF manually] で最適化を行う範囲を入力(3塩基単位)、または選択肢から指定
- ④ 3種類の条件を追加可能
  - Protect my cloning sites]:最適化時に変更したくない制限酵素サイトを指定可能
  - Protected Regions]:最適化時に変更したくない配列 領域を指定可能
  - [Motifs to Avoid]:最適化後の配列に含めたくない制
     限酵素サイトを指定可能(A)
- (B) [Fix Problems](問題点のみ最適化)または[Optimize](全長 を最適化)をクリックして最適化を実行、[Success]と表示され ればOK
- 6 複数配列を一括設定する場合は<u>P32</u>~を参照



**Thermo Fisher** 

SCIENTI

関連ページ:<u>P12, P13, P28, P32, P33</u>

### GeneArt Strings DNA Fragments Order カートに追加

- 価格・作業期間を確認して、注文ご希望のコンストラクト に☑
- ② [Download Summary]で配列情報を入手して確認
- ③ [Add to Cart] をクリックして注文手続きへと進む(詳細は [GeneArt人工遺伝子合成サービス ご注文マニュアル] のP8[6 カート内での価格・予想処理時間の確認]参照)

Optimize (optional)	Order		
Your Sequence Configuration	Your Products   Vertical Select all	Suggestions	JPY 133,028.00
			Project: 2022AADC3F Status: Draft
I HsActin_nuc	JPY 32,700 6 business days		Item Quantity DNA fragment 3 Cloned gene 1
ື່ HsActin_aa 1134 bp	Stitugs JPY 32,700 6 business days		Estimated production time 🟮 Strings - 6 business days Cloned Genes - 19 business days
២ nonString1 1054 bp	Shings JPY 32,700 6 business days		Email will be sent when Quality Assurance files are available in ThermoFisher Connect Kindly review the sequence(s) before ordering
ີງ nonString2 1054 bp	Ctoned genes JPY 34,928 19 business days	Fix/optimize to get string or Order as cloned gene	2 Download Summary 3 Add to Cart (4)

**Thermo Fisher** 

SCLEN

関連ページ:<u>P28</u>

# 複数遺伝子への一括設定適用・設定の保存 GeneArt Strings DNA Fragments編

#### 複数遺伝子への一括設定適用

Thermo Fisher SCIENTIFIC

 1番上のチェックボックスに☑を入れて、複数遺伝子を 一括選択

- 2 [Configure Selected] をクリック
- ③ 二通りの選択肢から選択
  - A) [Configure sequences manually]:新規に設定
  - B) [Apply project template]: 既存の雛形から選択
- (4) [Continue]をクリック
- 5 設定画面へ移動。次の2つのタブで各種設定が可能
- **(DNA Fragments Settings**]:
  - A) 5'/3'付加配列
- ⑦ [Optimization Settings]
  - A) 最適化生物種
  - B) 除外する配列

Add Sequence + Import & Save 4 Sequence(s) 2 Configure Selected Delete Selected Protein sequence(s) must have a host organism selected. You can do this for one or multiple sequences () 1 Host Organism 👙 Actions Status Sequence Name Туре Length 🕕 DNA/Protein Sequence 1 Delete 0 ofigure HsActin\_nuc DNA 1152 ATGGAATTCAAGCTTGGATCCTGTGAAGAAGAGGACAGCACTGCCTTGGTG. Cricetulus griseus Delete 0 MEFKLGSCEEEDSTALVCDNGSGLCKAGFAGDDAPRAVFPSIVGRPRHQGVM. HsActin\_aa PROTEIN 384 Homo sapiens 🔟 Delete 🖉 Co Ø HsActin\_nuc\_2 ATGGAATTCAAGCTTGGATCCTGTGAAGAAGAGGACAGCACTGCCTTGGTG. DNA 1152 Cricetulus griseus 🔟 Delete 🖉 Configure 0 HsActin\_aa\_2 PROTEIN 384 MEFKLGSCEEEDSTALVCDNGSGLCKAGFAGDDAPRAVFPSIVGRPRHQGVM Homo sapiens **Choose Configure Options** Х Configure Multiple Sequence(s) Х 5 Informational: The bulk configuration will only be applied to the selected sequences that are qualified. Kindly review the changes for individual sequence after saving Informational: Fields that cannot be configured in bulk are not displayed. Changes will **3**A Configure sequences manually be applied to the selected sequences. **③**B Apply project template 6 (7 4 Select the DNA Fragments Settings Optimization Settings Choose Configure Options Х 1 Informational: The bulk configuration will only be applied to the selected sequences that are qualified. Kindly review the changes for individual sequence after sav Configure sequences manually Apply project temp Select the project template (**3**)**B** Q search Cancel Continue Manual 230830

#### 関連ページ:<u>P12, P13, P27, P29</u>

#### Thermo Fisher

### 各種設定・雛形として保存

Optimization Settings	
3' Region(Optional) 🟮	(-
ТААТАА	
(or)	
& 3' region 🜖	
ete 3' Region	
	Iptimization Settings 3' Region(Optional) TAATAA (or) & 3' region ete 3' Region

① [5'/3' Region (Optional)]: 付加配列

	DNA Fragments Settings Optimiz	ation S	ettings						
	Host Organism for optimization 🔒				<b>^</b>				
기	Homo sapiens 👻								
)	Optimise the sequence upon Import & Save 0								
D	Motifs to Avoid Define motifs to avoid in your sequence. Selected sites will be excluded during optimization.								
	Standard Motifs Custom Motifs								
	Click on the motifs to add or remove 🕤								
	Standard Motifs		Standard Motifs to Avoid						
	Hind Q		Search motifs	Q	-				

- [Host Organism...]:最適化生物種を選択。最 適化しない場合は[No optimization]を選択
- ② [Optimize the sequence...]: ☑を入れると Import & Save 中に最適化を実施
- ③ [Motifs to Avoid]: 最適化配列から除外する配列

Hind	Q	Search motifs	Q
HindIII [AAGCTT]	*	BamHI [GGATCC]	*
		ECORI [GAATTC]	
	_		_
4		4	
		(or)	
Choose to delete any e	xisting motiffs to a	avoid 📵	
Delete motiffs to a	void		
)			
Save settings as proje	ect template 🖯		
roject Template Name *			
emplate Manual 230831			

- ① [Save settings as...]: ☑を入れると雛形として保存
- ② [Project Template Name]: 雛形の名前を入力\*1
- ③ [Save changes]: クリックして一括適用・雛形保存

\*1:全角文字・スペースはエラーの原因となりますので、ご注意ください。

#### 関連ページ: <u>P12</u>, <u>P13</u>, <u>P27</u>, <u>P29</u>

### 複数遺伝子への一括設定適用の完了

- (Import & Save)をクリック、Orderへ移動、この間に最 適化を実行(P28へ続く)

eque	nce(s)						Add Sequence	<ul> <li>Import &amp; Sav</li> </ul>
Con	figure Selected 🔞 Delete	Selected	Protein sequence	e(s) must have a	host organism sele	ected. You can do this for one or <b>multiple sequences ()</b>		
/	Actions	Status	Sequence Name	Туре 🍦	Length	DNA/Protein Sequence		Host Organism 🛓
	🖻 Delete 🖉 Configure	0	HsActin_nuc	DNA	1152	ATGGAATTCAAGCTTGGATCCTGTGAAGAAGAGGACAGCACTGC	CTTGGTG	Homo sapiens
/	🖻 Delete 🖉 Configure	0	HsActin_aa	PROTEIN	384	MEFKLGSCEEEDSTALVCDNGSGLCKAGFAGDDAPRAVFPSIVGRF	PRHQGVM	Homo sapiens
/	🕲 Delete 🖉 Configure	0	HsActin_nuc2	DNA	1152	ATGGAATTCAAGCTTGGATCCTGTGAAGAAGAGGACAGCACTGC	CTTGGTG	Homo sapiens
/	💼 Delete 🖉 Configure	0	HsActin_aa2	PROTEIN	384	MEFKLGSCEEEDSTALVCDNGSGLCKAGFAGDDAPRAVFPSIVGR	PRHQGVM	Homo sapiens

#### 関連ページ:<u>P28</u>

### ー括設定雛形の確認・修正

Thermo Fisher

- ① [Project Templates]から[DNA Fragments] を選択
- [Template Name]から該当の雛形をクリックして内容を 確認
- ③ 修正する場合は[Configure]または[Edit]をクリック
- ④ 新規に作成する場合は[Add New Template]をクリック



GeneArt	Project Templates - DNA F	Fragments
Welcome	Add or edit project templates with your the optimizati	ion settings to enable fast and easy submission of future projects.
A Home	Actions Template N	Name ≑ Last Modified ≑
Grders & Drafts	Delete 🖉 Configure	07-Sep-2023
₩ Project Templates ▲		
Cloned Genes		
DNA Fragments		
CopoArt	Back	
GeneAlt	test	3 Edit Template
Welcome		
<b>•</b>	DNA Fragments Settings Opt	imization Settings
Home	5' Region:	3' Region:
📜 Orders & Drafts	GCCACC	ТАДТАД
👭 Project Templates	; ▲	
Cloned Genes		
DNA Fragments		
GonoArt	•	
General	Edit Project Template	
Welcome	Save optimization settings as a project tem	plate that can be applied to future DNA fragments projects.
♠ Home	Project Template Name *	
	test	
Orders & Drafts	DNA Fragments Settings Optimiz	ation Settings
👫 Project Templates 🔺		
Cloned Genes	5' Region(Optional)	3' Region(Optional) 🕄
	GCCACC	TAGTAG

Thermo Fisher



# 過去に合成した遺伝子のプラスミド精製

36 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

#### Thermo Fisher

## 再注文 対象コンストラクトの選択

- た枠の対象となるプロジェクトの+をクリック(A)、プロ ジェクトとコンストラクトの名前・IDで検索も可能(B)
- ② 対象となるコンストラクトに☑(A)、配列情報をダウン ロードも可能(B)
- ③ [Add]をクリックして右枠へ追加
- ④ 対象となるコンストラクトに☑
- 5 [Import & Save]をクリック



37 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

Plasmid Reorder - Instant Designer	
Select the constructs for reordering	Cancel Import & Save
Constructs that are not available for reorder are displ	
Trojects and Constructs Search by project name or id Q	Constructs (Ottems) Search by construct name or id Q
Project Name 🝦 Project Id 💠 Last Updated ≑	Construct Name 🖕 Construct Id 🖕
+ 09/15/2022	No constructs selected
+ 09/15/2022	
<b>_</b>	
My Projects and Constructs Search by construct name or id Q	Constructs Otems Search by construct name or id Q
	✓ Construct Name
	No constructs selected
Construct Name   Construct Id   Action	10 20
▲         ▲	He frage
✓         ★         Remove	
Select the constructs for reordering	Cancel Import & Save
Constructs that are not available for reorder are displayed in grey 0	
My Projects and Constructs Search by construct name or id Q	ructs (211ems) Search by construct name or id Q
<ul> <li>⊀ Back to projects list</li> </ul>	Construct Name 🖕 Construct Id 🖕
2822AAKHOD 2022AAKHOD 09/15/2022	Internet and the second sec
Construct Name 🖕 Construct Id 🖕 Action	
× * *	
✓ End to the second	move

#### Thermo Fisher s c i e n t i f i c

## 再注文 プラスミド精製スケール・バッファーの選択

- ① ご希望のスケール・バッファーを選択(詳細は<u>P18</u>参照)
- 2 選択後[Confirm]をクリックして全てのコンストラクトに一括設定
- ③ 個別設定も可能
- ④ [Continue]をクリック



関連ページ:<u>P18</u>

#### 再注文 カートに追加

#### ① 価格・作業期間を確認

- [Download Summary]で配列情報を入手して確認
- ③ [Add to Cart] をクリックして注文手続きへと進む(詳細 は[GeneArt人工遺伝子合成サービス\_ご注文マニュア ル]のP8[6カート内での価格・予想処理時間の確認]参 照)

Plasmid Preparation Configuration	Your Product	JPY 74,200.00
		Project: 2022AAAAOR Status: Draft
TE, 1mg yield (megaprep), Transfection Grade	JPY 25300.00 6 business days	Item Quant Plasmid Preparation
3524 bp     TE. 1-10 mo vield (oloapreo). Transfection Grade	JPY 45900.00 6 business days	Estimated production time 0 Plasmid Preparation - 6 business days
		Email will be sent when Quality Assurance files are available in ThermoFisher Connect
		Kindly review the sequence(s) before ordering

関連ページ:

Thermo Fisher SCIENTIFIC

# プロジェクト概要の確認

40 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

### Drafts 設定中のプロジェクト

 [Drafts]で設定中・未注文のプロジェクトを表示可 能

② [Project Name]から各プロジェクト設定画面へ 移動可能

- ③ プロジェクトのタグも表示
- ④ 様々な項目でソート可能
- 5 プロジェクトのタグで抽出も可能
- 6 Project ID, Construct ID, Construct Nameなど で検索も可能

GeneArt	Drafts 214			Filter by tag	- Search	by project name or proj	ect id	Q
Home	G Add/Modify tag	G Remove tag		4				
In Florie	Action	Project Name \$	Project Id	# Sequences	Project Type \$	Last modified ÷	Status ¢	
<ul> <li>Orders &amp; Drafts</li> <li>Drafts</li> </ul>		Test_Strings_01	2024ABAL5F	2	DNA Fragments	14-Nov-2024	Draft	<b>^</b>
Open Orders Past Orders Offline Orders	📄 📋 Delete	Test01	2024ABHQ5D	2	Cloned Genes	14-Nov-2024	Draft	

#### 関連ページ:<u>P3</u>

#### Thermo Fisher

## **Open Orders** 進行中のプロジェクト

- [Open Orders]で進行中のプロジェクトを表示可能
- 2 [+]をクリックしてコンストラクト情報を表示 可能
- ③ 各項目でソート可能
- Actionの[<sup>1</sup>]をクリックすると次の3点が可能
  - A) 納品用配列情報ファイルのダウンロード
  - B) プロジェクトのOrderタブへの移動
  - c) Eメール通知の設定

GeneArt **Open Orders I** Q Search by project name or project id Filter by tag -3 Welcome Export all orders to Excel Order Est. Last A Home Project Name ± Project ID \$ Access 🚹 PO Number 🚖 Status 🚖 Action ٥ Update Shipment Number 2 📜 Orders & Drafts + St. Hannel Drafts Open Orders 🚹 +S damage . Past Orders 2 Concernance of the Institute of the Inst -09/22/2022 O In progress 10/04/2022 Construct ID Construct Name Product Type Est. Shipment Status 👙 10/04/2022 In progress .... Council per-10/04/2022 In progress -Donal gave COMPANY NAMES 10/04/2022 In progress ..... Correct pers 224.417162



関連ページ:<u>P3</u>

#### Past Orders 納品済みプロジェクト

- [Past Orders]で納品済みプロジェクトを表示可 能
- 2 [+]をクリックしてコンストラクト情報を表示可能
- ③ 各項目でソート可能
- ④ Actionの[<sup>1</sup>]をクリックすると次の2点が可能
  - A) 納品用配列情報ファイルのダウンロード
  - B) プロジェクトのOrderタブへの移動

	GeneArt <sup>Welcome</sup>	Past Orders 100		Filter by tag •	Search by project name or project id	Q
	A Home	Project Name   Project ID	Access 1 PO	Number    Order Number	er Status ≑	Action
	Drafts	+	Actorest		Pages 171000	1.1
1	Open Orders 1 Past Orders Offline Orders	+	A local	10000	Frager - 10.000	
2						
		2117482204	100.000	Shipped 1	1/13/2017 :	
	1 item(s) Construct Name	$\stackrel{\mathbb{A}}{_{\!$	Construct ID 🖕 St	atus 🛓		
		Crewit gene	00.07	Shipped		



#### 関連ページ:<u>P3</u>

#### 納品形態·出荷

- 人工遺伝子合成・サブクローニング:凍結乾燥DNA5µg
- プラスミド精製:DNA溶液(濃度1 mg/mL)
- GeneArt Strings DNA Fragments: 凍結乾燥DNA >200 ng
- 容器:チューブ(96ウェルプレート納品も可能、ご希望の場合は<u>専用納品依頼書にてご依頼</u>)
- 合成・構築が完了したコンストラクトから順に出荷。完了予定日前でも出荷。