

CAUTION

Before using this product, please read the Limited Use License statement below:

Important Limited Use License information for pFUSEN-Lucia-mG2AFc

The purchase of the pFUSEN-Lucia-mG2AFc vector conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes.

The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes.

Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic, or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research.

If the purchaser is unwilling to accept the limitations of this limited use statement, InvivoGen is willing to accept return of the product with a full refund. The product must be returned in resaleable condition. For information on purchasing a license to this product for purposes other than research, contact InvivoGen, 3950 Sorrento Valley Blvd. Suite 100, San Diego, CA 92121. Tel: 858-457-5873 Fax: 858-457-5843.

TECHNICAL SUPPORT

Toll free (US): 888-457-5873
Outside US: (+1) 858-457-5873
Europe: +33 562-71-69-39
E-mail: info@invivogen.com
Website: www.invivogen.com

 **InvivoGen**
3950 Sorrento Valley Blvd. Suite 100
San Diego, CA 92121 - USA

pFUSEN-Lucia-mG2AFc

Plasmid designed for Lucia::Fc fusion to the N-terminus of a protein of interest

Catalog # pfcn-lcmg2a

For research use only

Version 20K09-MM-v36

PRODUCT INFORMATION

Content:

- 20 μ g of pFU SEN-Lucia-mG2AFc plasmid provided as lyophilized DNA
- 1 ml of Zeocin™ (100 mg/ml)

Storage and Stability:

- Product is shipped at room temperature.
- Lyophilized DNA should be stored at -20°C and is stable 3 months.
- Resuspended DNA should be stored at -20°C and is stable up to 1 year.
- Store Zeocin™ at 4 °C or at -20 °C. The expiry date is specified on the product label.

Quality control:

- Plasmid construct has been confirmed by restriction analysis and sequencing.
- Plasmid DNA was purified by ion exchange chromatography and lyophilized.

GENERAL PRODUCT USE

pFUSEN-Fc / pFUSEN-Lucia-Fc is a family of plasmids developed to facilitate the construction of Fc-fusion proteins where the immunoglobulin G (IgG) Fc-domain is fused to the N-terminus of the protein of interest.

pFUSEN-Fc / pFUSEN-Lucia-Fc plasmids yield high levels of Fc-fusion proteins. The level of expression is usually in the μ g/mL range. They can be transfected in a variety of mammalian cells, including myeloma cell lines, CHO cells, monkey COS cells and human embryonic kidney (HEK) 293 cells, cells that are commonly used in protein purification systems.

pFUSEN-Lucia-mG2AFc plasmid allows the production of Lucia-Fc fusion proteins. This plasmid can be used to make recombinant Lucia-Fc fusion proteins or can be used as a transfection control in experiments with other pFUSEN-Fc constructs. Quantification of Lucia-Fc expression can be determined utilizing InvivoGen's QUANTI-Luc™ (rep-qlc1 or rep-qlc2).

A choice of cloning sites is provided to allow flexibility in the design of the fusion linker: either use pFUSEN linker, or bring forth your own linker with the protein of interest.

InvivoGen provides pFUSEN-Lucia-Fc vectors featuring Fc regions from different species and isotypes. In mouse, the IgG2a isotype is available. In humans, three options are available: IgG1, IgG1e2, or IgG2. The Fc region mediates effector functions, such as antibody-dependent cellular cytotoxicity (ADCC) and complement-dependent cytotoxicity (CDC). IgG isoforms exert different levels of effector functions. The engineered IgG1e2 contains mutations in the FcRn binding sites leading to higher FcRn binding affinity and reduced pH dependence.

PLASMID FEATURES

- **Lucia luciferase** is a secreted coelenterazine-utilizing luciferase reporter protein with advantageous characteristics when associated with Fc-fusion proteins. It possesses superior carrier ability for excellent secretion of the chimeric protein. It provides a simple means to screen for recombinant clones and it minimally affects the activity of the protein of interest.

- **Mouse IgG2a-Fc** : The Fc region comprises the CH2 and CH3 domains of the IgG2a heavy chain, with the hinge region. The last amino acid (lysine) of the Fc region has been replaced by an alanine for better fusion result. The Fc region of mouse IgG2a mediates high ADCC and CDC.

- **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1 α (EF-1 α) core promoter¹ and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat². The EF-1 α promoter exhibits a strong activity and yields long lasting expression of a transgene *in vivo*. The R-U5' has been coupled to the EF-1 α core promoter to enhance stability of RNA.

- **Cloning sites & fusion linker:** The protein of interest can be cloned either as a BamHI—NheI fragment, or as a BsiWI—NheI fragment. With BamHI cloning, the protein of interest will be separated from the Fc-domain by a flexible linker (Gly4Ser dimer).

With BsiWI cloning, the flexible linker will not be retained, allowing for a different fusion design. The provided cloning sites are compatible with many other enzymes, thus facilitating cloning.

- **SV40 pAn:** the Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA³.

- **ori:** a minimal *E. coli* origin of replication to limit vector size, but with the same activity as the longer Ori.

- **CMV enh / hFerL prom:** This composite promoter combines the human cytomegalovirus immediate-early gene 1 enhancer and the core promoter of the human ferritin light chain gene. This ubiquitous promoter drives the expression of the Zeocin™-resistance gene in mammalian cells.

- **EM2KC** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*. EM2KC is located within an intron and is spliced out in mammalian cells.

- **Zeo:** Resistance to Zeocin™ is conferred by the *Sh ble* gene from *Streptoalloteichus hindustanus*. The same resistance gene confers selection in both mammalian cells and *E. coli*.

- **β Glo pAn:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription⁴.

1. Kim DW *et al.* 1990. Use of the human elongation factor 1 alpha promoter as a versatile and efficient expression system. 91(2):217-23.

2. Takebe Y. *et al.* 1988. SR alpha promoter: an efficient and versatile mammalian cDNA

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3622-3480

E-mail: info@invivogen.com

expression system composed of the simian virus 40 early promoter and the R-U5 segment of human T-cell leukemia virus type 1 long terminal repeat. Mol Cell Biol. 8(1):466-72.

3. Carswell S. & Alwine JC. 1989. Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. Mol Cell Biol. 9(10):4248-58.

4. Yu J. & Russell JE. 2001. Structural and functional analysis of an mRNP complex that mediates the high stability of human beta-globin mRNA. Mol Cell Biol. 21(17):5879-88.

METHODS

Plasmid resuspension

Quickly spin the tube containing the lyophilized plasmid to pellet the DNA. To obtain a plasmid solution at 1 µg/µl, resuspend the DNA in 20 µl of sterile H₂O. Store resuspended plasmid at -20 °C.

Plasmid amplification and cloning

Plasmid amplification and cloning can be performed in *E. coli* GT116 or in other commonly used laboratory *E. coli* strains, such as DH5α.

Zeocin™ usage

This antibiotic can be used for *E. coli* at 25 µg/ml in liquid or solid media and at 50-200 µg/ml to select Zeocin™-resistant mammalian cells.

RELATED PRODUCTS

Product	Catalog Code
Zeocin™	ant-zn-1
QUANTI-Luc™	rep-qlc1

[TECHNICAL SUPPORT](#)

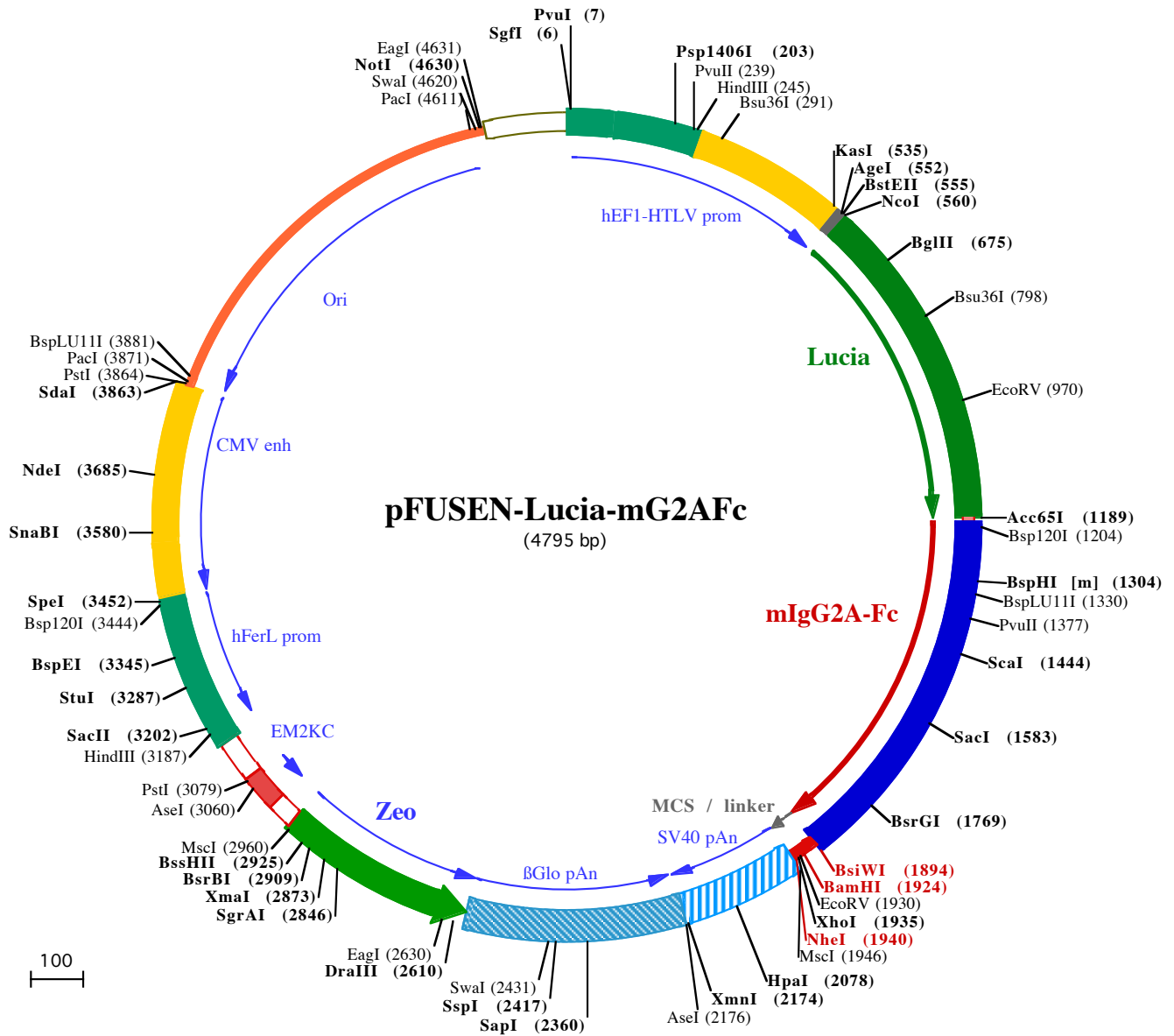
InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3622-3480

E-mail: info@invivogen.com



PvuI (7)
SgfI (6)
1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203) **HindIII (245)** **PvuII (239)** **Bsu36I (291)**
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACACGCCGCCGCCCTACCTGAGGCC
301 GCCATCCACGCCGGTTGAGTGCAGTTCGCCGCTCCCGCCTGTGGTGCCTCCTGAAGCTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC
401 GGGCCTTTGTCCGGCGCTCCCTTGAGGCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTCTCAACTCTACGTCTTTGTTTCGTTT

NcoI (560) **BstEII (555)**
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCGCTACCTGAGATCACCGGTACCATGAAATCAAGGTGCTGTTTGCCTCATCTGTATTGC
1 M E I K V L F A L I C I A
601 TGTGCTGAGGCAAACCCACTGAAATCAATGAAGACCTCAATATAGCTGCTGTGGCCTCCAACCTTGGCCACCACAGATCTTGAGACTGACCTGTTCCACC
13 V A E A K P T E I N E D L N I A A V A S N F A T T D L E T D L F T
701 AACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACTCCCCCAGATGTCC
47 N W E T M N V I S T D T E Q V N T D A D R G K L P G K K L P P D V
801 TGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTGGTGCACAAGAGGCTGCCTCATTGGCTCTCCACATTAAGTGACCCCTAAGATGAAGAAATTTAT
80 L R E L E A N A R R A G C T R G C L I C L S H I K C T P K M K K F I
901 CCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCCAGAGATTCTGGCTTCAAGGAT
113 P G R C H T Y E G E K E S A Q G G I G E A I V D I P E I P G F K D
1001 AAGGAGCCACTGGACCACTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGCTGTCTGAAGGGCCTTGCCAAATGTCCAGTCTCTGACCTCC
147 K E P L D Q F I A Q V D L C A D C T T G C L K G L A N V Q C S D L
1101 TGAAGAAGTGGCTTCCCCAGAGGTGACCACTTTTCCAGCAAGATTGAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGACAGAGGTACCCAGGCC
180 L K K W L P Q R C T T F A S K I Q G R V D K I K G L A G D R G T E P
Bsp120I (1204)
1201 CAGAGGGCCACAATCAAGCCCTGTCTCCATGCAAAATGCCAGCACCTAACCTTTGGTGGACCATCCGTCTTCATCTTCCCTCCAAGATCAAGGAT
2 R G P T I K P C P P C K C P A P N L L G G P S V F I F P P K I K D
BspHI [m] (1304) **BspLU11I (1330)** **PvuII (1377)**
1301 GTACTCATGATCTCCCTGAGCCCATAGTCACATGTGTGGTGGATGTGAGCGAGGATGACCCAGATGTCCAGATCAGCTGGTTTGTGAACAACGTGG
36 V L M I S L S P I V T C V V V D V S E D D P D V Q I S W F V N N V
1401 AAGTACACACAGCTCAGACACAAACCCATAGAGAGGATTACAACAGTACTCTCCGGTGGTCACTGCCCCTCCCATCCAGCACCAGGACTGGATGAGTGG
69 E V H T A Q T Q T H R E D Y N S T L R V V S A L P I Q H Q D W M S G
ScaI (1444) **SacI (1583)**
1501 CAAGGAGTCAAATGCAAGGTCAACAACAAGACCTCCAGCGCCATCGAGAGAACCATCTCAAACCAAAGGGTCAAGAGCTCCACAGGTATAT
102 K E F K C K V N N K D L P A P I E R T I S K P K G S V R A P Q V Y
1601 GTCTTGCCTCCACCAGAAGAAGAGATGACTAAGAAACAGGTCACTCTGACCTGCATGGTCACAGACTTCATGCCTGAAGACATTTACGTGGAGTGGACCA
136 V L P P P E E E M T K K Q V T L T C M V T D F M P E D I Y V E W T
BsrGI (1769)
1701 ACAACGGGAAAACAGAGCTAAACTACAAGAACACTGAACAGTCTGGACTCTGATGGTCTTACTTTCATGTACAGCAAGCTGAGAGTGGAAAAGAAGAA
169 N N G K T E L N Y K N T E P V L D S D G S Y F M Y S K L R V E K K N
BsiWI (1894)
1801 CTGGTGGAAAGAAATAGCTACTCTGTTCACTGTTGAGTGGTCCAGGGTCTGCACAATCACCACACGACTAAGAGCTTCTCCGGACTCCGGGTGACAGTACG
202 W V E R N S Y S C S V V H E G L H N H H T T K S F S R T P G A
1 R T
EcoRV (1930) **NheI (1940)**
1901 GGTGGTGGCGGTAGCGGTGGTGGCGGATCCGATATCTCGAGCTAGCTGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATG
3 G G G G S G G G G S D I S S •
HpaI (2078)
2001 CAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAACAAGTTAAACAACAACATTGCATTC
AseI (2176) **XmnI (2174)**
2101 ATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGTGTGTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTAATTCTAAAATACAGCATAGC
2201 AAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTT
SapI (2360)
2301 GCAGCCTCACCTTCTTTTCATGGAGTTTAAAGATATAGTGTATTTTCCCAAGTTTGAAGTACTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCCA

2401 **SspI (2417)** **SwaI (2431)**
 CATTCCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCTT

2501 CATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTC
 125 D Q

2601 **DraIII (2610)** **EagI (2630)**
 GCTCCTCTGCCACAAAGTGCACGCAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTCATGGCCGGCCCGGA
 122 E E A V F H V C N G A P D R L A F E R G W P Q E G I E T M A P G S

2701 GCGTCCCGGAAGTTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCCGCGCACCCACACCCAGGCCAGGGTGTGTCCGGCACCACC
 89 A D R F N T S V V E S W E A Y L E D L G R V W V W A L T N D P V V

2801 **SgrAI (2846)** **XmaI (2873)**
 TGGTCTGGACCGCGCTGATGAACAGGGTACGTCGTCGCCGACACACCGGCGAAAGTCTCCTCCACGAAGTCCCGGAGAACCCGAGCCGGTTCGGTCC
 55 Q D Q V A S I F L T V D D R V V G A F D D E V F D R S F G L R D T W

2901 **BsrBI (2909)** **BssHIII (2925)** **MscI (2960)**
 AGAACTCGACCGCTCCGGCGACGTCGCGCGGGTGAGCACCAGGAAACGGCAGTGGTCAACTTGGCCATGATGGCTCTCctgtcaggagaggaagagaag
 22 F E V A G A V D R A T L V P V A S T L K A M

3001 aaggttagtacaattgCTATAGTGAGTTGATTATACTATGCAGATATACTATGCCAATGATTAATTGTCAAACCTAGGGCTGCagggttcattagtccac
 AseI (3060) PstI (3079)

3101 ttttctgcactgccccatctctgcccaccctttccaggcatagacagtcagtgacttacCAAACCTACAGGAGGGAGAAGGCAGAAGCTTGAGACAG
 HindIII (3187)

3201 **SacII (3202)** **StuI (3287)**
 ACCCGCGGACCGCCGAACCTGCGAGGGGACGTGGCTAGGGCGGCTTCTTTTATGTTGCGCGGCCCTCGAGGCAGGGCGCTCGGGGAGCCCTAGCGGCC

3301 **BspEI (3345)**
 AATCTGCGGTGGCAGGAGCGGGGCCGAAGCCGTGCCTGACCAATCCGGAGCACATAGGAGTCTCAGCCCCCGCCCCAAAGCAAGGGGAAGTCACGCG

3401 **SpeI (3452)** **BspI20I (3444)**
 CCTGTAGCGCCAGCGTGTGTGAAATGGGGCTTGGGGGGTTGGGGCCCTGACTAGTCAAAACAAACTCCATTGACGTCAATGGGGTGAGACTTGGGA
 SpeI (3452)

3501 AATCCCCGTGAGTCAAACCGCTATCCACGCCCATTTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAG
 SnaBI (3580)

3601 TAGGAAAGTCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTG
 NdeI (3685)

3701 ATGTACTGCCAAGTGGGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCATTATTGAC

3801 **PacI (3871)** **PstI (3864)** **SdaI (3863)** **BspLU11I (3881)**
 GTCAATGGGCGGGGTCGTTGGGCGGTACGCCAGGCGGGCCATTTACCGTAAGTTATGTAACCGCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCA
 SdaI (3863)

3901 GCAAAAGGCCAGAACCGTAAAAAGCCGCTTGTGCGCTTTTCCATAGGCTCCGCCCTGACGAGCATCAAAAAATCGACGCTCAAGTCAGAGG

4001 TGGCGAAACCCGACAGGACTATAAAGATACCAGGCTTTCCCTGGAAGTCCCTCGTGGCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGT

4101 CCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAGCTGGGCTGTGTGCA

4201 CGAACCCCCGTTACGCCGACCGCTGCGCCTTATCCGTAACATATCGTCTTGTAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACT

4301 GGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCT

4401 GCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGAAGCA

4501 GCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAAACGAAAACCTCACGTTAAGGGATTTG

4601 **EagI (4631)** **PacI (4611)** **SwaI (4620)** **NotI (4630)**
 GTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTAATCGTAACTAAC

4701 ATACGCTCTCCATCAAAACAAAACGAAACAAAACAAACTAGCAAAATAGGCTGTCCCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA