

CAUTION

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

Important Limited Use License information for pFUSEN-Lucia-hG1Fc

The purchase of the pFUSEN-Lucia-hG1Fc 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, 10515 Vista Sorrento Parkway San Diego, CA 92121 USA. Tel: 858-457-5873 Fax: 858-457-5843.

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



pFUSEN-Lucia-hG1Fc

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

Catalog # pfcn-lchg1

For research use only

Version 20K09-MM-v36

PRODUCT INFORMATION

Content:

- 20 μ g of pFUSEN-Lucia-hG1Fc 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-hG1Fc 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 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.
- **Human IgG1-Fc** : The Fc region comprises the CH2 and CH3 domains of the IgG1 heavy chain, with the hinge region. The first cysteine of the hinge has been replaced by a serine to prevent detrimental disulfide bridges. The last amino acid (lysine) of the Fc region has been replaced by an alanine for better fusion result. Human IgG1 displays high ADCC and CDC, and is the most suitable for therapeutic use against pathogens and cancer cells.
- **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*.
- **βGlo pAn**: The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription⁴.

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



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 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

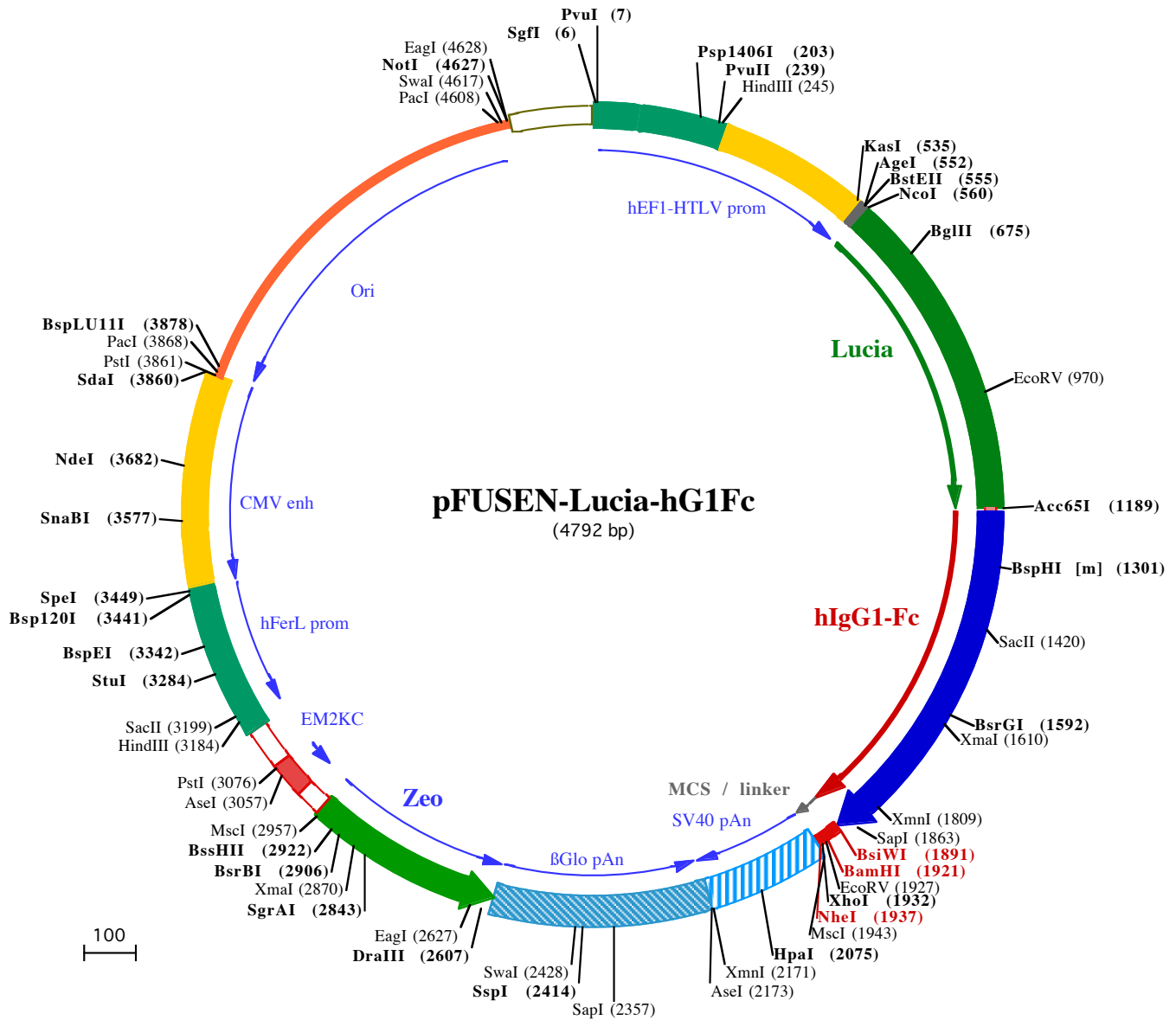
InvivoGen USA (Toll-Free): 888-457-5873

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 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTACTGGCTCCGCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)
Psp1406I (203) **PvuII (239)**
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACCGCGCCGCCGCCCTACCTGAGGCC
301 GCCATCCACGCCGGTTGAGTGCAGTCTGCCGCTCCCGCCTGTGGTGCCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC
401 GGGCCTTTGTCCGGCGCTCCCTTGGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGTCTTTGTTTCGTTT

NcoI (560) **BstEII (555)**
KasI (535) **AgeI (552)**
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTACCATTGAAATCAAGGTGCTGTTTGCCTCATCTGTATTGC
1 M E I K V L F A L I C I A
BglIII (675)
601 TGTGCTGAGGCAAACCCACTGAAATCAATGAAGACCTAATATAGCTGCTGTGGCCTCCAACCTTGGCCACCACAGATCTTGAGACTGACCTGTTCCACC
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 TGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTGGTGCACAAGAGGCTGCCTCATTTGCCTCTCCACATTAAGTGCACCCTAAGATGAAGAAATTTAT
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
EcoRV (970)
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 AAGGAGCCACTGGACCAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGGCCTTGCCAATGTCCAGTGTCTGACCTCC
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
Acc65I (1189)
1101 TGAAGAAGTGGCTTCCCCAGAGGTGTACCACTTTTGCAGCAAGATTCAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGAGGTACCGAGCC
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
hinge Cys changed to Ser (1207)
1201 CAAATCTAGTGACAAAACCTCACACATGCCACCCTGCCAGCACCTGAACTCCTGGGGGACCGTCAGTCTTCTCTTCCCCCAAAACCAAGGACACC
2 K S S D K T H T C P P C P A P E L L G G P S V F L F P P K P K D T
BspHI [m] (1301)
1301 CTCATGATCTCCCGACCCCTGAGGTCACATGCGTGGTGGACGTGAGCCACGAAGACCCTGAGGTCAAGTCAACTGGTACGTGGACGGCGTGGAGG
36 L M I S R T P E V T C V V V D V S H E D P E V K F N W Y V D G V E
SacII (1420)
1401 TGCATAATGCCAAGACAAAGCCGCGGGAGGAGCAGTACAACAGCACGTACCGTGTGGTCAAGCTCCTACCGTCTGCACCAGGACTGGCTGAATGGCAA
69 V H N A K T K P R E E Q Y N S T Y R V V S V L T V L H Q D W L N G K
BsrGI (1592)
1501 GGAGTACAAGTGAAGGTCTCCAACAAAGCCCTCCAGCCCCATCGAGAAAACCATCTCAAAGCCAAAGGGCAGCCCCGAGAACCAGGTGTACACC
102 E Y K C K V S N K A L P A P I E K T I S K A K G Q P R E P Q V Y T
XmaI (1610)
1601 CTGCCCATCCCGGAGGAGATGACCAAGAACCAGGTGACGCTGACCTGCTGGTCAAAGGCTTCTATCCAGCGACATCGCCGTGGAGTGGGAGAGCA
136 L P P S R E E M T K N Q V S L T C L V K G F Y P S D I A V E W E S
1701 ATGGGAGCCGGAGAACAACACTACAAGACCACGCTCCCGTGTGGACTCCGACGGCTCCTTCTTCTTACAGCAAGCTCACCGTGGACAAGAGCAGGTG
169 N G Q P E N N Y K T T P P V L D S D G S F F L Y S K L T V D K S R W
BsiWI (1891)
XmnI (1809) **SapI (1863)** **Lys changed to Ala (1888)**
1801 GCAGCAGGGGAAAGCTCTTCTCATGCTCCGTGATGCACGAGGCTCTGCACAACCACTACACGCAGAAGGCCTCTCCCTGTCTCCGGGTGCAAGTACGGGT
202 Q Q G N V F S C S V M H E A L H N H Y T Q K S L S L S P G A
1 R T G
EcoRV (1927) **NheI (1937)**
BamHI (1921) **XhoI (1932)** **MscI (1943)**
1901 GGTGGCGGTAGCGGTGGTGGCGGATCCGATATCTCGAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAG
4 G G G S G G G S D I S S •
HpaI (2075)
2001 TGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAAGTTAAACAACAATTCATT
AseI (2173)
XmnI (2171)
2101 TTATGTTTCAGGTTCAAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTAATTCTAAAATACAGCATAGCAA
2201 ACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTTTGCCAATGTGCATTAGCTGTTTGA
SapI (2357)
2301 GCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCCAAGGTTTGAAGTACTGCTCTTCAATTTCTTTATGTTTTAAATGCACTGACCTCCACAT

SspI (2414) SwaI (2428)
2401 TCCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCAT
2501 AATATCCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCCTGCT
125 • D Q E
DraIII (2607) EagI (2627)
2601 CCTCTGCCACAAAGTGCACGCAGTTGCCGGCCGGTCCGCGAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTCATGGCCGGCCCGGAGGC
121 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 A
2701 GTCCCGAAAGTTCGTGGACACGACCTCCGACCCTCGGCGTACAGCTCGTCCAGGCCGCGCACCCACACCCAGGCCAGGGTGTGTCCGGCACCACTGG
88 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 Q
SgrAI (2843) XmaI (2870)
2801 TCCTGGACCGCGCTGATGAACAGGGTCACGTCGTCGCCGACCACACGGCGAAGTCTCTCCACGAAGTCCCGGGGAGAACCAGCCGGTCCGGTCCAGA
54 D Q V A S I F L T V D D R V V G A F D E V F D R S F G L R D T W F
BsrBI (2906) BssHII (2922) MscI (2957)
2901 ACTCGACCGCTCCGGCGACGTCGCGCGGGTGAGCACCGGAACGGCACTGGTCAACTGGCCATGATGGCTCCTctgtcaggagaggaagagagaagaag
21 E V A G A V D R A T L V P V A S T L K A M
AseI (3057) PstI (3076)
3001 gttagtacaattgCTATAGTGAGTTGTATTATACTATGCAGATATACTATGCCAATGATTAATTGCAAACTAGGGCTGCAgggttcatagtgcacttt
HindIII (3184) SacII (3184)
3101 tcctgcaactgcccattctcctgcccacccttccaggcatagacagtcaagtgttacCAAACCTCACAGGAGGGAGAAGGCAGAAAGCTTTGAGACAGACC
StuI (3284)
3201 CGCGGGACCGCCGAACCTGCGAGGGGACGTGGCTAGGGCGGCTCTTTTATGGTGCGCCGCCCTCGGAGGCAGGGCGCTCGGGAGGCCCTAGCGGCCAAT
BspEI (3342)
3301 CTGCGGTGGCAGGAGGCGGGGCCGAAGCCGTGCTGACCAATCCGGAGCACATAGGAGTCTCAGCCCCCGCCCCAAAGCAAGGGGAAGTCACGCGCCT
SpeI (3449)
Bsp120I (3441)
3401 GTAGCGCCAGCGTGTGTGAAATGGGGCTTGGGGGGTGGGGCCCTGACTAGTCAAAACAAACTCCCATTGACGTCAATGGGGTGGAGACTTGAAAT
SnaBI (3577)
3501 CCCCCTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAG
NdeI (3682)
3601 GAAAGTCCATAAGGTCACTGACTGGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGATG
3701 TACTGCCAAGTGGGCAAGTTTACCCTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTC
PacI (3868)
PstI (3861)
SdaI (3860) BspLU11I (3878)
3801 AATGGGCGGGGTCGTTGGGCGGTCAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCCCTGCAGGTTAATTAAGAACATGTGAGCAAAAGCCAGCA
3901 AAAGGCCAGGAACCGTAAAAAGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGG
4001 CGAAACCCGACAGGACTATAAGATAACCAGGCGTTTTCCCTGGAAAGCTCCCTCGTGCCTCTCTGTTCGACCCTGCCGTTACCGGATACCTGTCCG
4101 CCTTTCTCCCTTCGGAAGCGTGGCGTTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGACGA
4201 ACCCCCCGTTCCAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGT
4301 AACAGGATTAGCAGAGCGAGGTATGTAGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCG
4401 CTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTTGTTTGAAGCAGCA
4501 GATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGGAAACGAAACTCACGTTAAGGGATTTTGGTC
EagI (4628)
PacI (4608) SwaI (4617) NotI (4627)
4601 ATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATA
4701 CGCTCTCCATCAAAACAAACGAAACAAACAAACTAGCAAAATAGGCTGTCCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA