

# pSELECT-zeo-LucSh

A plasmid encoding a synthetic CpG-free firefly Luc-Zeocin resistance fusion gene

Catalog code: psetz-lucsh

For research use only

Version 20K30-MM

## PRODUCT INFORMATION

### Content:

- 20 µg of pSELECT-zeo-LucSh 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 resuspended upon receipt and stored at -20°C. Lyophilized DNA is stable for 3 months at -20°C. Resuspended DNA is stable more than one year at -20°C.

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

pSelect-zeo plasmids contain genes that have been chemically synthesized. The DNA sequence of these genes was modified by optimizing the codon usage, reducing or eliminating the CpG motifs and avoiding secondary DNA structures without changing the amino acid sequence of the wild type proteins.

pSelect-zeo plasmids may be used:

**To subclone the synthetic gene into another vector.** To facilitate subcloning, the LucSh gene is flanked by two unique restriction sites: Nco I at the 5' end that encompasses the Start codon, and Nhe I at the 3' end.

**As a gene reporter plasmid.** pSelect-zeo is a mammalian expression plasmid selectable in *E. coli* and mammalian cells with Zeocin™, as the *Sh ble* gene in the second expression cassette is driven by the eukaryote CMV enhancer/promoter in tandem with the bacterial EM7 promoter.

## PLASMID FEATURES

### First expression cassette

• **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1alpha (EF-1α) core promoter<sup>1</sup> and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat<sup>2</sup>. The EF-1α promoter exhibits a strong activity and yields long lasting expression of a transgene *in vivo*. The R-U5' has been coupled to the EF-1α core promoter to enhance stability of RNA.

• **LucSh:** Synthetic LucSh fusion gene (LucSh-ΔCpG): InvivoGen has engineered a fusion between the firefly luciferase gene and the *Sh ble* gene conferring Zeocin™ resistance. Both genes have been modified and contain no CpG, whereas their wildtype counterparts contain 95 and 50 CpG motifs respectively. This fusion exhibits a higher luciferase activity and enables a better and faster selection of Zeocin™ resistant clones.

• **SV40 pAn:** the Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA<sup>3</sup>.

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

### Second expression cassette

• **CMV enh/prom:** The human cytomegalovirus immediate-early gene 1 promoter/enhancer was originally isolated from the Towne strain and was found to be stronger than any other viral promoters.

• **EM7** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*.

• **Zeo:** Resistance to Zeocin™ is conferred by the *Sh ble* gene from *Streptoalloteichus hindustanus*. The *Sh ble* gene is driven by the CMV enhancer/promoter in tandem with the bacterial EM7 promoter allowing 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 transcription<sup>4</sup>.

1. Kim, D.W. *et al.* (1990). *Gene* 2: 217-223.

2. Takebe, Y. *et al.* (1988). *Mol. Cell Biol.* 1: 466-472.

3. Carswell, S. & Alwine, J.C. (1989). *Mol. Cell Biol.* 10: 4248-4258.

4. Yu J & Russell JE. (2001). *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 H<sub>2</sub>O. Store resuspended plasmid at -20 °C.

### **Plasmid amplification and cloning**

Plasmid amplification and cloning can be performed in *E. coli* GT116 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..

## TECHNICAL SUPPORT

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

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

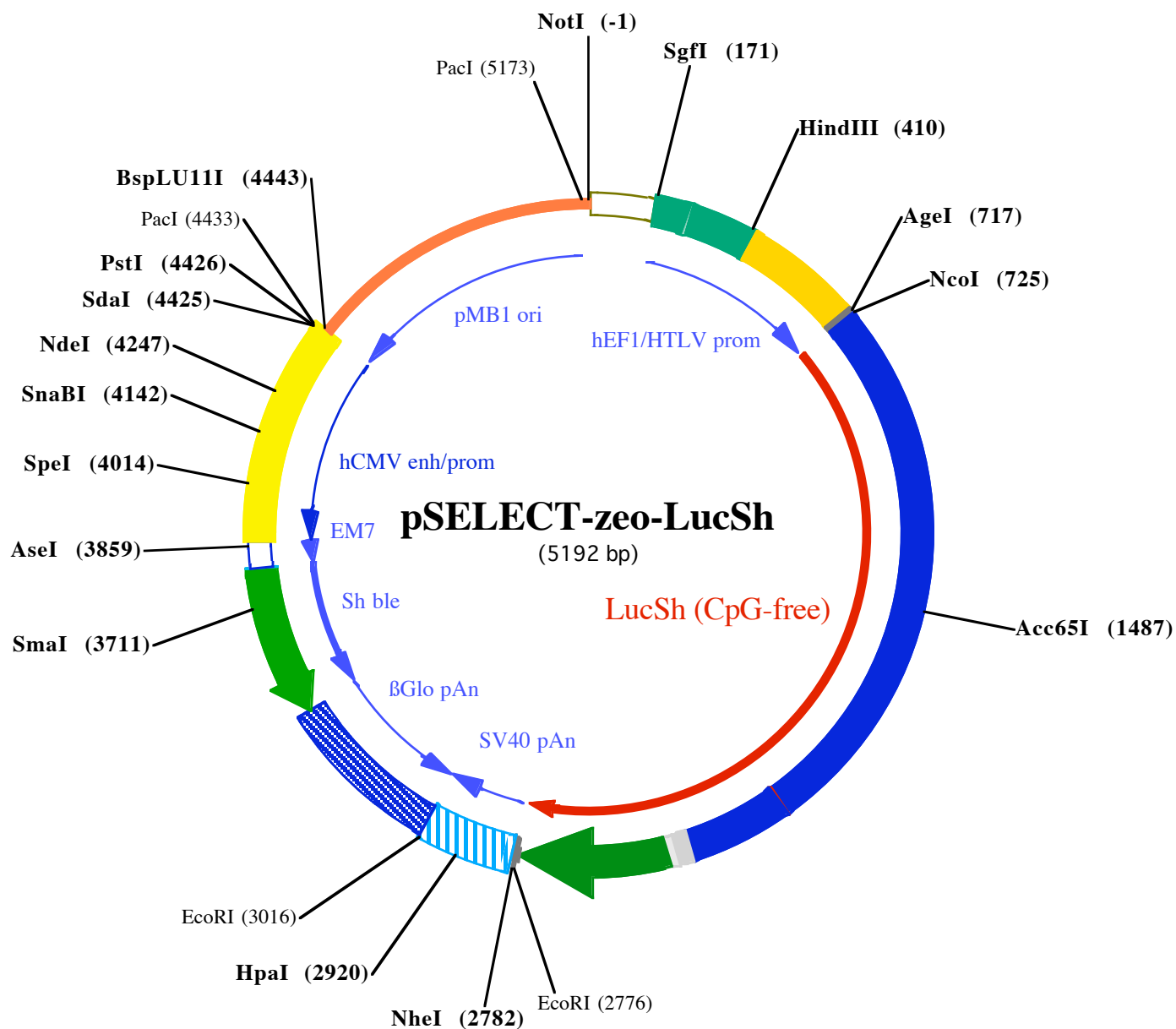
InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)



[www.invivogen.com](http://www.invivogen.com)



**NotI (-1)**

1 GCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACT  
68 AACATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAAATAGGCTGTCCCCAGTGCA

**SgfI (171)**

135 AGTGCAGGTGCCAGAACATTTCTCTATCGAAGGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGCAG  
202 AGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGGAGGGGTGCGCAATTGAACGGGTGCCTAGAG  
269 AAGGTGGCGCGGGTAAACTGGGAAAGTGATGTCGTGACTGGCTCCGCCTTTTTCCCGAGGGTGGG  
336 GGAGAACCGTATATAAGTGCAGTAGTCGCCGTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAA

**HindIII (410)**

403 CACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTCACGCGCCCGCCGCTACCTGAGGCCGCCA  
470 TCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAACTGCGTCCGCCGCTC  
537 TAGGTAAGTTTAAAGCTCAGGTCGAGACCGGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGAC  
604 TCAGCCGGCTCTCCACGCTTTGCCTGACCCTGCTTGCTCAACTCTACGTCTTTGTTTCGTTTTCTGT

**NcoI (725)**

**AgeI (717)**

671 TCTGCGCCGTTACAGATCCAAGCTGTGACCGGCGCCTACCTGAGATCACCGGTACCATGGAGGATG  
738 CCAAGAATATTAAGAAAGGCCCTGCCCATCTACCTCTGGAAGATGGCACTGCTGGTGAGCAACT  
4▶A K N I K K G P A P F Y P L E D G T A G E Q L  
805 GCACAAGGCCATGAAGAGGTATGCCCTGGTCCCTGGCACCATTGCCTTCACTGATGCTCACATTGAG  
26▶H K A M K R Y A L V P G T I A F T D A H I E  
872 GTGGACATCACCTATGCTGAATACTTTGAGATGTCTGTGAGGCTGGCAGAAGCCATGAAAAGATATG  
49▶V D I T Y A E Y F E M S V R L A E A M K R Y  
939 GACTGAACACCAACCACAGGATTGTGGTGTGCTCTGAGAACTCTCTCCAGTTCTTCATGCCTGTGTT  
71▶G L N T N H R I V V C S E N S L Q F F M P V L  
1006 AGGAGCCCTGTTCAATTGGAGTGGCTGTGGCCCCTGCCAATGACATCTACAATGAGAGAGAGCTCCTG  
93▶G A L F I G V A V A P A N D I Y N E R E L L  
1073 AACAGCATGGGCATCAGCCAGCCAAGTGTGGTCTTTGTGAGCAAGAAGGGCCTGCAAAAGATCCTGA  
116▶N S M G I S Q P T V V F V S K K G L Q K I L  
1140 ATGTGCAGAAGAAGCTGCCCATCATCCAGAAGATCATCATCATGGACAGCAAGACTGACTACCAGGG  
138▶N V Q K K L P I I Q K I I I M D S K T D Y Q G  
1207 CTTCCAGAGCATGTATACCTTTGTGACCAGCCACTTACCCCTGGCTTCAATGAGTATGACTTTGTG  
160▶F Q S M Y T F V T S H L P P G F N E Y D F V  
1274 CCTGAGAGCTTTGACAGGGACAAGACCATTGCTCTGATTATGAACAGCTCTGGCTCCACTGGACTGC  
183▶P E S F D R D K T I A L I M N S S G S T G L  
1341 CCAAAGGTGTGGCTCTGCCCCACAGAAGTGTGTGAGATTGAGCCATGCCAGAGACCCCATCTT  
205▶P K G V A L P H R T A C V R F S H A R D P I F  
1408 TGGCAACCAGATCATCCCTGACACTGCCATCCTGTCTGTGGTTCCATTCCATCATGGCTTTGGCATG  
227▶G N Q I I P D T A I L S V V P F H H G F G M

**Acc65I (1487)**

1475 TTCACAACACTGGGGTACCTGATCTGTGGCTTCAGAGTGGTGCTGATGTATAGGTTTGAGGAGGAGC  
250▶F T T L G Y L I C G F R V V L M Y R F E E E  
1542 TGTTTCTGAGGAGCCTACAAGACTACAAGATCCAGTCTGCCCTGCTGGTGCCCACTCTGTTTCAGCTT  
272▶L F L R S L Q D Y K I Q S A L L V P T L F S F  
1609 CTTTGCCAAGAGCACCTCATTGACAAGTATGACCTGAGCAACCTGCATGAGATTGCCTCTGGAGGA  
294▶F A K S T L I D K Y D L S N L H E I A S G G  
1676 GCACCCCTGAGCAAGGAGGTGGGTGAGGCTGTGGCAAAGAGGTTCCATCTCCAGGAATCAGACAGG  
317▶A P L S K E V G E A V A K R F H L P G I R Q  
1743 GCTATGGCCTGACTGAGACCACCTCTGCCATCCTCATCACCCCTGAAGGAGATGACAAGCCTGGTGC  
339▶G Y G L T E T T S A I L I T P E G D D K P G A  
1810 TGTGGGCAAGGTGGTTCCCTTTTTTGAGGCCAAGGTGGTGGACCTGGACACTGGCAAGACCCTGGGA  
361▶V G K V V P F F E A K V V D L D T G K T L G  
1877 GTGAACCAGAGGGGTGAGCTGTGTGTGAGGGGTCCCATGATCATGTCTGGCTATGTGAACAACCCTG  
384▶V N Q R G E L C V R G P M I M S G Y V N N P

1944 AGGCCACCAATGCCCTGATTGACAAGGATGGCTGGCTGCACTCTGGTGACATTGCCTACTGGGATGA  
406▶ E A T N A L I D K D G W L H S G D I A Y W D E  
2011 GGATGAGCACTTTTTTCATTGTGGACAGGCTGAAGAGCCTCATCAAGTACAAAGGCTACCAAGTGGA  
428▶ D E H F F I V D R L K S L I K Y K G Y Q V A  
2078 CCTGCTGAGCTAGAGAGCATCCTGCTCCAGCACCCCAACATCTTTGATGCTGGTGTGGCTGGCCTGC  
451▶ P A E L E S I L L Q H P N I F D A G V A G L  
2145 CTGATGATGATGCTGGAGAGCTGCCTGCTGCTGTTGTGGTTCTGGAGCATGGAAGACCATGACTGA  
473▶ P D D D A G E L P A A V V V L E H G K T M T E  
2212 GAAGGAGATTGTGGACTATGTGGCCAGTCAGGTGACCACTGCCAAGAAGCTGAGGGGAGGTGTGGTG  
495▶ K E I V D Y V A S Q V T T A K K L R G G V V  
2279 TTTGTGGATGAGGTGCCAAAGGGTCTGACTGGCAAGCTGGATGCCAGAAAGATCAGAGAGATCCTGA  
518▶ F V D E V P K G L T G K L D A R K I R E I L  
2346 TCAAGGCCAAGAAGGGTGGCAAACAATTGATCTCTGGAGCCAATGGAGTATGGCCAAGTTGACCAG  
540▶ I K A K K G G K Q L I S G A N G V M A K L T S  
2413 TGCTGTCCAGTGCTCACAGCCAGGGATGTGGCTGGAGCTGTTGAGTTCTGGACTGACAGGTTGGGG  
562▶ A V P V L T A R D V A G A V E F W T D R L G  
2480 TTCTCCAGAGATTTTGTGGAGGATGACTTTGCAGGTGTGGTCAGAGATGATGTCACCCTGTTTCATCT  
585▶ F S R D F V E D D F A G V V R D D V T L F I  
2547 CAGCAGTCCAGGACCAGGTGGTGCCTGACAACACCCTGGCTTGGGTGTGGGTGAGAGGACTGGATGA  
607▶ S A V Q D Q V V P D N T L A W V W V R G L D E  
2614 GCTGTATGCTGAGTGGAGTGGTGTCTCCACCAACTTCAGGGATGCCAGTGGCCCTGCCATGACA  
629▶ L Y A E W S E V V S T N F R D A S G P A M T  
2681 GAGATTGGAGAGCAGCCCTGGGGGAGAGAGTTTGCCTGAGAGACCCAGCAGGCAACTGTGTGCACT  
652▶ E I G E Q P W G R E F A L R D P A G N C V H

**NheI (2782)**

EcoRI (2776)

2748 TTGTGGCAGAGGAGCAGGACTGAGGATAAGAATTGCTAGCTGGCCAGACATGATAAGATACATTGA  
674▶ F V A E E Q D •  
2815 TGAGTTTGGACAAACCACAACACTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCT

**HpaI (2920)**

2882 ATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCATTTTA  
2949 TGTTTCAGGTTTCAGGGGGAGGTGTGGGAGGTTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTAT

EcoRI (3016)

3016 GGAATTCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCT  
▶ ◀  
3083 GAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACC  
3150 TTCTTTCATGGAGTTTAAGATATAGTGTATTTTCCCAAGGTTTGAAGTACTAGCTCTTCATTTCTTTATG  
3217 TTTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCA  
3284 TTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCC  
3351 CAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGC  
3418 TTCTAGCTTATCCTCAGTCCTGCTCCTCTGCCACAAAGTGCACGCAGTTGCCGGCCGGGTTCGCGCAG  
125▶ • D Q E E A V F H V C N G A P D R L  
3485 GCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTCATGGCCGGCCCGGAGGCGTCCCGGAAG  
107▶ A F E R G W P Q E G I E T M A P G S A D R F  
3552 TTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCCGCGCACCCACACCCAGGCCA  
84▶ N T S V V E S W E A Y L E D L G R V W V W A L  
3619 GGGTGTGTCCGGCACCACCTGGTCTGGACCGCGCTGATGAACAGGGTCACGTCGTTCCCGGACCAC  
62▶ T N D P V V Q D Q V A S I F L T V D D R V V

**SmaI (3711)**

3686 ACCGGCGAAGTCGTCTCCACGAAGTCCCGGAGAACCAGCCGGTCCGAGCCGGTCCAGAACTCGACCGCT  
40▶ G A F D D E V F D R S F G L R D T W F E V A  
3753 CCGGCGACGTCGCGCGCGGTGAGCACCGGAACGGCACTGGTCAACTTGGCCATGATGGCCCTCTAT  
17▶ G A V D R A T L V P V A S T L K A M ◀

**AseI (3859)**

3820 AGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGC

3887 GTCTCCAGCTTATCTGACGGTTCCTAAACGAGCTCTGCTTATATAGACCTCCCACCGTACACGCCT

**SpeI (4014)**

3954 ACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAG

4020 TCAAAACAAACTCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTA

**SnaBI (4142)**

4087 CCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGT

4154 ACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGCATAATGCCAGGCGGGCCATTTACCGTC

**NdeI (4247)**

4221 ATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGCAGTTTAC

4288 CGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCA

4355 TTATTGACGTCAATGGGGCGGGGTCGTTGGGCGGTCAGCCAGGCGGGCCATTTACCGTAAGTTATGT

**PacI (4433)**

**PstI (4426)**

**SdaI (4425)**

**BspLU11I (4443)**

4422 AACGCCTGCAGGTTAA TTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAA

4487 GGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCA

4554 AGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCCCTGGAAGCTCCCTCG

4621 TGCGCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCCTTTCTCCCTTCGGGAAGCGT

4688 GGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTTCGCTCCAAGCTGGGC

4755 TGTGTGCACGAACCCCCGTTCCAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCA

4822 ACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTA

4889 TGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTT

4956 GGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAAC

5023 AAACCACCGCTGGTAGCGGTGGTTTTTTTTGTTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATC

5090 TCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGG

**PacI (5173)**

5157 ATTTTGGTCATGGCTAGTTAATTAACATTTAAATC A