

STOP

Before opening this package, please read the Limited Use License statement below:

Important Limited Use License information for pCpGfree-vitroHLacZ

The purchase of the pCpGfree-vitroHLacZ 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 A, 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 A
San Diego, CA 92121 - USA

pCpGfree-vitroHLacZ

A LacZ expression plasmid completely devoid of CpG dinucleotides, selectable with hygromycin

Catalog # pcpvyth-lz

For research use only

Version # 09E11-MM

PRODUCT INFORMATION

Content:

- 20 µg of pCpGfree-vitroHLacZ plasmid provided as lyophilized DNA
- *E. coli* GT115 strain provided lyophilized on a paper disk
- 4 pouches of Fast-Media® Hygro (2 TB and 2 Agar)

Storage and Stability:

- Product is shipped at room temperature.
- Lyophilized DNA is stable for 12 months when stored at -20°C.
- Resuspended DNA is stable for 6 months when stored at -20°C.
- Bacteria should be stored at -20°C and are stable up to 1 year.
- Store Fast-Media® Hygro at room temperature. Fast-Media® pouches are stable 18 months when stored properly.

Quality control:

- Plasmid construct has been confirmed by restriction analysis and sequencing.
- Plasmid DNA was purified by ion exchange chromatography and lyophilized.
- Viability of the lyophilized bacteria upon resuspension has been verified.

GENERAL PRODUCT USE

pCpGfree-vitro plasmids represent innovative tools to study the effects of CpG dinucleotides in a number of applications. DNA vaccination exploits the immunostimulatory character of certain CpG motifs to prime and boost the immune response. However, these immunostimulatory CpG motifs are antagonized by CpG dinucleotides in certain distinct base contexts, termed neutralizing CpG motifs. Both types of CpG motifs are usually present in plasmidic DNAs, and therefore may lead to an unfavorable immune response. pCpGfree-vitro is the ideal tool to overcome this problem, and may be used to study the effects of these two types of CpG motifs by adding them in different configurations to the pCpGvitro backbone.

CpG dinucleotides are key elements in a number of cellular functions associated with chromatin. Many recent findings have altered our vision of chromatin and its role in the regulation of cellular processes such as transcription regulation, DNA replication and repair, cell cycle control, and cell aging. Several large multisubunit complexes, consisting of methyl-CpG binding (MBD) proteins and histone deacetylases, have been implicated in the regulation of chromatin dynamics. These complexes are recruited to methylated CpG dinucleotides by DNA methyl transferases (DNMTs) and induce chromatin remodelling. However the specific roles of these complexes are still to be explored. Due to the absence of CpG dinucleotides within its backbone, pCpGfree-vitro is not the target of DNMTs and thus MBD proteins. Therefore, it provides a useful model to study the other proteins involved in these complexes, in particular the histone deacetylases. It can also be used to analyze the effects of CpG methylation on the regulation and duration of gene expression.

PLASMID FEATURES

pCpGfree-vitro is a new family of expression vectors completely devoid of CpG dinucleotides that are selectable in mammalian cells. Similarly to the other pCpGfree plasmids (i.e. pCpGfree-lacZ, pCpGfree-mcs, and pCpGfree-siRNA), all the elements required for replication and selection of the plasmids in bacteria, and gene expression in mammalian cells have been modified to remove all CpG dinucleotides.

- **Composite CpG-free promoter** combining the mouse CMV enhancer, the human elongation factor 1 α core promoter and 5'UTR containing a synthetic intron (I 126). This composite promoter yields high and ubiquitous expression of the LacZ gene.
- **LacZ** encodes β -galactosidase an enzyme that catalyzes the hydrolysis of X-Gal, producing a blue precipitate that can be easily visualized under a microscope. This CpG-free allele of the *lacZ* reporter gene can be easily subcloned and replaced by a gene of interest.

- **CpG-free polyadenylation signals (pAn):** The polyadenylation signals utilized are CpG-free versions of the SV40 late and human β -globin polyadenylation signals. These polyA enable efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA.
- **CpG-free matrix attached regions (MARs)** are AT-rich sequences that are able to form barriers between independent expression cassettes.
- **CpG-free Hygromycin resistance gene (hph- Δ CpG):** The CpG-free Hygro^R gene is active both in *E. coli* and mammalian cells.
- **CpG-free SV40 promoter** works in tandem with a bacterial promoter located within a synthetic intron (I-EC2K). This composite promoter drives the expression of the resistance gene in both mammalian cells and *E. coli*.
- **CpG-free *E. coli* R6K gamma origin of replication:** This origin is activated by the R6K specific initiator protein π , encoded by the *pir* gene. Expression of the *pir* gene is necessary for the replication and amplification of pCpGvitro plasmids. *E. coli* GT115 strain expresses a *pir* mutant gene that allows higher plasmid copy number.

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.

Reconstitution of *E. coli* GT115 strain

Use sterile conditions to do the following:

- 1- Reconstitute *E. coli* GT115 by adding 1 ml of LB medium in the tube containing the paper disk. Let sit for 15 minutes. Mix gently by inverting the tube several times. Let sit 5 more minutes.
- 2- Streak bacteria taken from this suspension on a LB agar plate.
- 3- Place the plate in an incubator at 37°C overnight.
- 4- Isolate a single colony and grow the bacteria in *E. coli* growth medium.
- 5- Prepare competent cells utilizing your preferred protocol.

Preparation of Fast-Media Hygro

Fast-Media® is a **fast and convenient** way to prepare liquid and solid media for bacterial culture by using only a microwave. Fast-Media® Hygro can be ordered separately [#fas-hg-1 (TB), #fas-hg-s (Agar)].

Method:

- 1- Pour the contents of a Fast-Media® pouch into a clean borosilicate glass bottle or flask.
- 2- Add 200 ml of distilled water to the flask
- 3- Heat in a microwave on MEDIUM power setting (about 400Watts), until bubbles start appearing (approximately 3 minutes). **Do not heat a closed container. Do not autoclave Fast-Media®.**
- 4- Swirl gently to mix the preparation. **Be careful, the bottle and media are hot, use heatproof pads or gloves and care when handling.**
- 5- Reheat the media for 30 seconds and gently swirl again. Repeat as necessary to completely dissolve the powder into solution. But be careful to avoid overboiling and volume loss.
- 6- Let agar medium cool to 45°C before pouring plates. Let liquid media cool to 37°C before seeding bacteria.

Note: Do not reheat solidified Fast-Media® as the antibiotic will be permanently destroyed by the procedure.

References:

1. Wu F. *et al.* 1995., A DNA segment conferring stable maintenance on R6K gamma-origin core replicons. *J Bacteriol.* 177(22):6338-45.
2. Bode J. *et al.*, 1996. Scaffold/matrix-attached regions: topological switches with multiple regulatory functions. *Crit Rev Eukaryot Gene Expr.* 6(2-3):115-38.

TECHNICAL SUPPORT

Toll free (US): 888-457-5873

Outside US: (+1) 858-457-5873

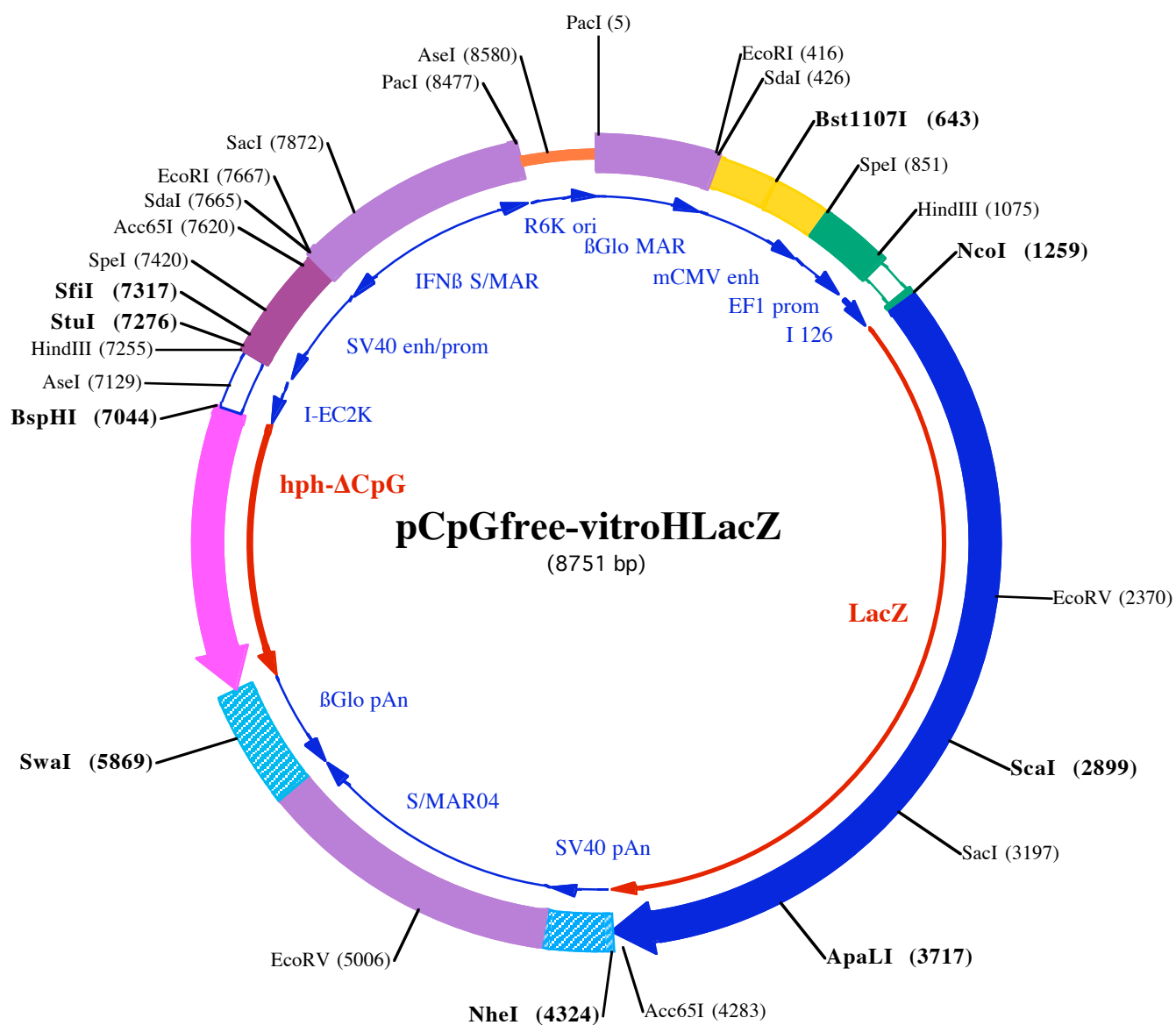
Europe: +33 562-71-69-39

E-mail: info@invivogen.com

Website: www.invivogen.com



3950 Sorrento Valley Blvd. Suite A
San Diego, CA 92121 - USA



225

PacI (5)
 1 TTAATTAATTAATCTCTAAGGCATGTGAAGTGGCTGTCTTGGTTTTTCATCTGTACTTCATCTGCTACCTCTGTGACCTGAAACATATTTATAATTCCAT
 101 TAAGCTGTGCATATGATAGATTTATCATATGTATTTTCTTAAAGGATTTTGTGAAGAACTAATTGAATTGATACCTGTAAAGTCTTTATCACACTACCC
 201 AATAAATAATAATCTCTTTGTCAGCTCTCTGTTCTATAAATGTACAAGTTTTATTGTTTTAGTGGTAGTGATTTATTCTCTTTCTATATATAT
 301 ACACACACATGTGTGCATTCAATAATATATACAATTTTTATGAATAAAAAATTATTAGCAATCAATATTGAAAACCACTGATTTTGTGTTATGTGAGCAA

 SdaI (426)
 EcoRI (416)
 401 ACAGCAGATTAAAGGAATTCCTGCAGGAGTCAATGGGAAAAACCATTGGAGCCAAGTACACTGACTCAATAGGGACTTTCCATTGGGTTTTGCCAGT
 501 ACATAAGGTCAATAGGGGTGAGTCAACAGGAAAGTCCATTGGAGCCAAGTACATTGAGTCAATAGGGACTTTCCAATGGGTTTTGCCAGTACATAAG

 Bst1107I (643)
 601 GTCAATGGGAGGTAAGCCAATGGGTTTTTCCATTACTGACATGTATACTGAGTCAATTAGGGACTTTCCAATGGGTTTTGCCAGTACATAAGGTCAATA
 701 GGGGTGAATCAACAGGAAAGTCCATTGGAGCCAAGTACACTGAGTCAATAGGGACTTTCCATTGGGTTTTGCCAGTACAAAAGGTCAATAGGGGGTGA

 SpeI (851)
 801 GTCAATGGGTTTTTCCATTATTGGCACATACATAAGGTCAATAGGGGTGACTAGTGGAGAAGAGCATGCTTGAGGGCTGAGTGCCCTCAGTGGGCAGA
 901 GAGCACATGCCCCACAGTCCCTGAGAAGTTGGGGGAGGGGTGGGCAATTGAACCTGGTGCCTAGAGAAGGTGGGGCTTGGGTAACTGGGAAAGTGATGT

 HindIII (1075)
 1001 GGTGTAAGTCCACCTTTTCCCCAGGGTGGGGGAGAACCATATATAAGTGCAGTAGTCTCTGTGAACATTCAAGCTTCTGCCTTCTCCCTCCTGTGA
 1101 GTTTGtaagtcactgactgtctatgcctgggaaggggtgggcaggagatggggcagtgaggaaaagtggcactatgaacceTGCAGCCCTAGCTAGAc

 NcoI (1259)
 1201 aattgtactaaccttcttctcttctcctcctgcagGTTGGGTACAGTAGCTTCCACCATGGACCCTGTTGTGCTGCAAAGGAGAGACTGGGAGAACC
 1301 CTGGAGTGACCCAGCTCAACAGACTGGCTGCCACCCTCCCTTTGCCTCTTGGAGGAACCTCTGAGGAAGCCAGGACAGACAGGCCAGCCAGCAGCTCAG
 1401 GTCTCTCAATGGAGAGTGGAGGTTTGCCTGGTTCCTGCCCTGAAGCTGTGCCTGAGTCTTGGCTGGAGTGTGACCTCCAGAGGCTGACACTGTTGTG
 1501 GTGCCAGCAACTGGCAGATGCATGGCTATGATGCCCCATCTACACCAATGTCACTACCCATCACTGTGAACCCCTTTTGTGCCACTGAGAACC
 1601 CCACTGGCTGCTACAGCTGACCTTCAATGTTGATGAGAGCTGGCTGCAAGAAGGCCAGACCAGGATCATCTTTGATGGAGTCAACTCTGCCTTCCACCT
 1701 CTGGTGCAATGGCAGGTGGGTTGGCTATGGCCAAGACAGCAGGCTGCCCTCTGAGTTTGACCTCTCTGCCTTCTCAGAGCTGGAGAGAAGCGTGGCT
 1801 GTCATGGTGCTCAGGTGGTCTGATGGCAGCTACCTGGAAGACCAAGACATGTGGAGGATGTCTGGCATCTTCAGGGATGTGAGCCTGCTGCACAAGCCCA
 1901 CCACCCAGATTTCTGACTTCCATGTTGCCACCAGGTTCAATGATGACTTCAGCAGAGCTGTGCTGGAGGCTGAGGTGCAGATGTGTGGAGAACTCAGAGA
 2001 CTACCTGAGAGTCACAGTGAGCCTCTGGCAAGGTGAGACCCAGGTGGCTCTGGCACAGCCCCCTTTGGAGGAGAGATCATTGATGAGAGAGGAGGCTAT
 2101 GCTGACAGAGTCACCTGAGGCTCAATGTGGAGAACCCCAAGCTGTGGTCTGCTGAGATCCCCAACCTCTACAGGGCTGTTGTGGAGCTGCACACTGCTG
 2201 ATGGCACCTGATTGAAGCTGAAGCCTGTGATGTTGGATTGAGAGAAGTCAGGATTGAGAATGGCCTGCTGCTGCTCAATGGCAAGCCTCTGCTCATCAG

 EcoRV (2370)
 2301 GGGAGTCAACAGGCATGAGCACCACCTCTGCATGGACAAGTGATGGATGAACAGACAATGGTGCAAGATATCTGCTAATGAAGCAGAACAACTTCAAT
 2401 GCTGTCAAGTGCTCTCACTACCCCAACCACCCTCTCTGGTACACCCTGTGTGACAGGTATGGCCTGTATGTTGTTGATGAAGCCAACATTGAGACACATG
 2501 GCATGGTGCCCATGAACAGGCTCACAGATGACCCAGGTGGCTGCCTGCCATGTCTGAGAGAGTGACCAGGATGGTGCAGAGAGACAGGAACACCCCTC
 2601 TGTGATCATCTGGTCTCTGGGCAATGAGTCTGGACATGGAGCAACCATGATGCTCTCTACAGGTGGATCAAGTCTGTTGACCCAGCAGACCTGTGCAG
 2701 TATGAAGGAGGTGGAGCAGACACCACAGCCACAGACATCATCTGCCCATGTATGCCAGGGTTGATGAGGACCAGCCCTTCCCTGCTGTGCCAAGTGA

 ScaI (2899)
 2801 GCATCAAGAAGTGGCTCTCTCTGCCTGGAGAGACCAGACCTCTGATCCTGTGTGAATATGCACATGCAATGGGCAACTCTCTGGGAGGCTTTGCCAAGTA
 2901 CTGGCAAGCCTTCAGACAGTACCCAGGCTGCAAGGAGGATTTGTGTGGGACTGGGTGGACCAATCTCTCATCAAGTATGATGAGAATGGCAACCCTGG
 3001 TCTGCCTATGGAGGAGACTTTGGTGACACCCCAATGACAGGCAGTTCTGCATGAATGGCCTGGTCTTTGACAGAGGACCCCTCACCTGCCCTCACAG

 SacI (3197)
 3101 AGGCCAAGCACCAGCAACAGTTCTTCCAGTTCAGGCTGTCTGGACAGACCATTGAGGTGACATCTGAGTACCTCTTCAGGCACTCTGACAATGAGTCTCT
 3201 GCACTGGATGGTGGCCCTGGATGGCAAGCCTCTGGCTTCTGGTGAAGTGCCTCTGGATGTGGCCCTCAAGGAAAGCAGCTGATTGAACTGCCTGAGCTG
 3301 CCTCAGCCAGAGTCTGCTGGACAACGTGGCTAACAGTGAGGGTGGTTCAGCCCAATGCAACAGCTTGGTCTGAGGCAGGCCACATCTCTGCATGGCAGC

3401 AGTGGAGGCTGGCTGAGAACCTCTCTGTGACCGCTGCTGCCTCTCATGCCATCCCTCACCTGACAACATCTGAAATGGACTTCTGCATTGAGCTGGG

3501 CAACAAGAGATGGCAGTTCAACAGGCAGTCTGGCTTCTGTCTCAGATGTGGATTGGAGACAAGAAGCAGCTCCTACCCCTCTCAGGGACCAATTACCC

3601 AGGGCTCCTCTGGACAATGACATTGGAGTGTCTGAGGCCACCAGGATTGACCCAAATGCTTGGGTGGAGAGGTGGAAGGCTGCTGGACACTACCAGGCTG

ApaLI (3717)

3701 AGGCTGCCCTGCTCCAGTGCACAGCAGACACCCCTGGCTGATGCTGTTCTGATCACCACAGCCCATGCTTGGCAGCACCAAGGCAAGACCCCTGTTTCATCAG

3801 CAGAAAGACCTACAGGATTGATGGCTCTGGACAGATGGCAATCACAGTGGATGTGGAGGTTGCCTCTGACACACCTCACCTGCAAGGATTGGCCTGAAC

3901 TGTCAACTGGCACAGGTGGCTGAGAGGGTGAACCTGGCTGGGCTTAGGCCCTCAGGAGAACTACCCTGACAGGCTGACAGCTGCCTGCTTTGACAGGTGGG

4001 ACCTGCCTCTGTCTGACATGTACACCCCTTATGTGTTCCCTTCTGAGAATGGCTGAGGTGTGGCACCAGGAGCTGAACTATGGTCCTACCAGTGGAG

4101 GGGAGACTTCCAGTTCAACATCTCCAGGTACTCTCAGCAACAGCTCATGAAACCTCTCAGGCACCTGCTCCATGCAGAGGAGGGAACCTGGCTGAAC

Acc65I (4283)

4201 ATTGATGGCTCCACATGGGCATTGGAGGAGATGACTCTTGGTCTCCTTCTGTGTCTGCTGAGTTCAGTTATCTGCTGGCAGGTACCACTATCAGCTGG

NheI (4324)

4301 TGTGGTGCCAGAAGTAAACCTGAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTT

4401 TATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACCAAGTTAACAAACAATTGCATTCATTTTATGTTTCAGGTT

4501 CAGGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTGAGGCCCACTGTGTTTCATCTTACAGATGGAATACTG

4601 ACATTGAGAGGAGTTAGTTAACTTGCTAGGTGATTGAGCTAATAAGTGAAGAAAGATTTCATCAAGGTGATTGATTCTGAAGCCTGTGCTAATCA

4701 CATTACCAAGCTACAACCTTCAATTATAATAATAAGTCAGCTTTCAAGGGCCTTTCAGGTGCTCTGCACCTTCTACAAGCTGTGCCATTAGTGAACAC

4801 AAAATGAGCCTTCTGATGAAGTAGTCTTTTCATTATTTAGATATTAGAACAATAAAATCTTAGCTGCCAGCTGATTGAAGGCTGGGACAAAATTCAAA

4901 CATGCATCTACAACAATATATATCTCAATGTTAGTCTCCAAATCTATTGACTTCAACTCAAGAGAATATAAAGAGCTAGTCTTTATACACTCTTTAAGG

EcoRV (5006)

5001 TATGATATCATCTGGAAAGTAACAAAATTGATGCAAATTTGAATGAACTTTATCATGGTGATTTACACAATGTGTTTCTTCCCTGCAATGTATTTCT

5101 TTCTCTAATTCCTTCCATTTGATCTTTCATACACAATCTGGTTCTGATGATGTTTTTGGATGCACCTTTCAACTCCAAAAGACAGAGCTAGTTACTTT

5201 CTTCTGGTGCTCCAAGCACTGTATTTGTATCTGTATTCAAGCCCTTGAATATTGTACTGGATCATTATTTACCTCTAGGATGGCTTCCCAGGCAA

5301 CTTGTGTTCCACCAGAGACTACATTTGTATCTTGTTGACCTTTGAACTTCCACCAGTGTCTAAAAATAATATGTATGCAAAATTACTTGCTATGAGAAT

5401 GTATAATTAACAATATAAAAAGGAGAAGCAAGGAGAGAAACACAGGTGTGATTTGTGTTTGTGTGCTTAAAGGCAGTGTGAAAAGGAAGAAATGCC

5501 ATTTATAGTGAGGAGACAAAGTTATATTACCTCTTATCTGGCTTTAAGGAGATTTGCTGAGCTAAAAATCCTATATTCATAGAAAAGCCTTACCTGAG

5601 TTGCCAATACCTCAATTCTAAAATACAGCATAGCAAACTTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGC

5701 ATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGACGCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCCAAGGTTGAACTAGCTCTT

SwaI (5869)

5801 CATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAAATATTAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTT

5901 TTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCAATATCCCCAGTTTAGTAGTTGACCTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAG

6001 CAAGAAAGCTCTAGCTAATATTCACTTCCTGGCTCTGGGCTTGTGGAGGGTCTCCTGTTTCCAGAGTCAGCCAGAAGTTCAACACATCCATCAGTCCAA

342 ●●● Gl uLysAl aArgProArgThr Ser ProArgArgAsnGl ySerAspAl aLeuVal Gl uVal CysGl yAspThr TrpV

6101 ACAGCAGCAGACCTTCTTGCAATTTGAGTTCTTCCAACAGTTCCAGCACCAGACCTCACAATGGCATCACATCTTCTTGTGCCCAAGCAGCATCATCAA

315 aAl aAl aSerArgArgAl aI l eGl nThrArgGl yVal Thr Gl yAl aGl ySerArgVal I l eAl aAspCysArgGl yGl nAl aTrpAl aAl aAspAspPh

6201 AGTTTCCATCAACCAGAGATTGATACAGTTGGTCCAGGCCAATTCTGAGCATGTAGGCTCTCAGTCTGGGGGAACCAAGCAGGTTCTGGGTGCTCTTCTTC

282 eAsnGl yAspVal LeuSer Gl nTyrLeuGl nAspLeuGl yI l eArgLeuMetTyrAl aArgLeuArgProSer Gl yAl aLeuGl uProHis sArgArgGl u

6301 AAAATATCTTGTGTTGTTCCATGCAAGCCAGCCAAGGCTCTCAAAAAAAATGTTGGCAACCTCATATTGAGAATCTCCAACATGGCTTCAGACCAAG

249 ePheTyrArgThr Gl nGl nGl uMetCysAl aLeuTrpProArgTrpPhePheI l eAsnAl aVal Gl uTyrGl nSerAspGl yPheMetAl aGl uSerTrpA

6401 TCAATGACTGCGATTCTGCCATTGCTGTGCGAAGCATTGTTGCTTCCAAAATCAGCATGGACCAGGTGCTGACTTCAGGACAATCTTCTGCCCAAC

215 spl l eValAl aThr I l eArgGl yAsnAspThr LeuVal AsnAsnSer Gl yPheAspAl aHisVal l eLeuHis sArgVal Gl uProCysAspGl uAl aTrpLe

6501 GCATGAGTTTCATCCAGTGCTGAGCAACAGAAAGCAGAACTGTGTCTCATCCATCAGCTCTGCCAGTGATAGACATGAGGATCAGCAATGGCACAATGAA

182 uMetLeuGl uAspLeuAl aGl nAl aVal l eSerAl aSerVal ThrAspAspMetVal l eThrGl nTrpHis sTyrVal l eHis sProAspAl aI l eAl aCysI l ePhe

6601 ATCCCTCCAAGTGGTGACTGACCAATGCCCTGGGAGCCAAAGGACCAATCCAGAGGTTTGGCTCAGATCAGCTGCTGCAATGGCATCCATTCCTTCA

149 AspArgTrpThr Thr TyrGl nGl yI l eGl yGl nProGl yPheProGl yPheGl ySer Thr Gl nSerLeuAspAl aAl aAl aI l eAl aAspMetAl aGl uA

6701 GCAACAGGTTGCAGAACAGCTGGCAGCTCAGTTTCAGGGAGATCCTGGAGAGTGACTCTTGTGCTCTTCTGCTGATGAGTAGGTTGAGAGATTCAGAAA

115 l aValProGl nLeuValAl aGl nLeuGl uThr Gl uProLeuAspGl nLeuThrVal Gl yGl nAl aArgArgSer I l eCysTyrThrLeuSerGl uSerPh

6801 ATTCTCAATGTCCAGAATCTTGAATTTGGCAGAGCAGCAGAGGCAAGTGTCTGTAACATATCTGTCTTTGTAAAAACCATCAGCACAAGAAATTGAC

82 eGl uGl yI l eAspLeuVal Gl uProI l eProLeuAl aAl aSerAl aPheHis sArgTyrVal l eTyrArgAspLysTyrPheGl yAspAl aCysSerAsnVal

6901 CCTCAGAACATAACCTCTTCTTCCACATCAAAAGAAAAGGCTCTGCTTCTTCCACCTTCAGACAGCTGCATGAGATCAGAAACAGAAATCAAAATTTTCA

49 ArgLeuVal l eTyrGl yArgGl yGl yVal AspPheSer PheAl aArgSer Gl uGl uGl yGl uSerLeuGl nMetLeuAspSerVal l eSerAspPheLysGl uI

7001 **ATGAGAACTTCTCAACAGAAGTTGCTGTCAAGTTTCAGTTTCAT**^{BspHI (7044)}**GATGGCTCCTCctgtcaggagaggaaagagaagaaggttagtacaattgCTAT**
 15 **↓** **I**eLeu**P**he**L**ys**G**lu**V**al**S**er**T**hr**A**l**a**Thr**L**eu**G**lu**P**ro**L**ys**L**ys**M**et

7101 **AGTGAGTTGTATTATACTATGCTTATGATTAATTGTCAAAGCTAGGGCTGC**^{AseI (7129)}**Agggttcatagtgccacttttctgcactgccccatctcctgccaccct**

7201 **ttcccaggcatagacagtcagtgacttacCAAAC**^{HindIII (7255)}**CACAGGAGGAGAAGGCAG**^{StuI (7276)}**AAGCTTTTTC****CAAAAGCCTAGGCCTCCAAAAAGCCTCCTCACTAC**

7301 **TTCTGGAATAGCTCAGAGGCCcAGGgGGCCTgGGCCTCTGCATAAATAAAAAAATTAGTCAGCCTGGGGctggggtgggggcaggggtggggggccaac**^{SfiI (7317)}

7401 **tgggCAGGGGTGGGGGGCCACTAGTGGGACTATGGTTGCTGACTAATTGAGATGCATGCTTTGCATACTTCTGCCTGCTGGGAGCCTGGGGACTTTCCA**^{SpeI (7420)}

7501 **CACCTGGTTGCTGACTAATTGAGATGCATGCTTTGCATACTTCTGCCTGCTGGGAGCCTGGGACTTTCCACACCCTAACTGACACACATTCCACAGCT**

7601 **GGTTCCTTCAGCCTCAGAGGTACCTAACCAAGTTCCTCTTTCAGAGGTTATTTTCAGGCCCTGCAGGAATTCAGTCAATATGTTACCCCCAAAAAGCTG**^{Acc65I (7620)}^{EcoRI (7667)}^{SdaI (7665)}

7701 **TTTGTTAACTTGTC AACCTCATTCTAAAATGTATATAGAAGCCAAAGACAATAACAAAAATATTCTTGTAACAAAATGGGAAAGAATGTTCCACTA**

7801 **AATATCAAGATTTAGAGCAAAGCATGAGATGTGTGGGATAGACAGTGAGGCTGATAAAATAGAGTAGAGCTCAGAAACAGACCCATTGATATATGTAAG**^{SacI (7872)}

7901 **TGACCTATGAAAAAATATGGCATTTTACAATGGGAAATGATGGTCTTTTTCTTTTTTAGAAAAACAGGGAAATATATTTATATGTA AAAAATAAAAGG**

8001 **GAACCCATATGTCATACCATAACACAAAAAAATTCAGTGAATTATAAGTCTAAATGGAGAAGGCAAAACCTTTAAATCTTTAGAAAAATAATATAGAAG**

8101 **CATGCCATCAAGACTTCAGTGTAGAGAAAAATTTCTTATGACTCAAAGTCCTAACCAAAAGAAAAGATTGTTAATTAGATTGCATGAATATTAAGACTT**

8201 **ATTTTTAAAAATTA AAAAACCTTAAGAAAAGTCAGGCCATAGAATGACAGAAAAATTTTGCAACACCCAGTAAAGAGAATTGTAATATGCAGATTATAA**

8301 **AAAGAAGTCTTACAAATCAGTAAAAAATAAACTAGACAAAAATTTGAACAGATGAAAGAGAAACTCTAAATAATCATTACACATGAGAAACTCAATCTC**

8401 **AGAAATCAGAGAACTATCATTGCATATACACTAAATTAGAGAAATATTTAAAGGCTAAGTAACATCTGTGGCTTAATTAAGTTATCCTAGGAAACCTTAA**^{PaeI (8477)}

8501 **AACCTTTAAAGCCTTATATATTCTTTTTTTCTTATAAAACTTAAACCTTAGAGGCTATTTAAGTTGCTGATTATATTAATTTTATTGTTCAAACAT**^{AseI (8580)}

8601 **GAGAGCTTAGTACATGAACATGAGAGCTTAGTACATTAGCCATGAGAGCTTAGTACATTAGCCATGAGGGTTTAGTTCATTAACATGAGAGCTTAGTA**

8701 **CATTAACATGAGAGCTTAGTACATACTATCAACAGGTTGAAGTCTGCTGATT**