

STOP

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

Important Limited Use License information for pVITRO2-blasti-Lucia/SEAP

The purchase of the pVITRO2-blasti-Lucia/SEAP 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 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



pVITRO2-blasti-Lucia/SEAP

A multigenic plasmid for high levels of expression of the Lucia luciferase and SEAP reporter genes

Catalog code: pvitro2-blucsp

<https://www.invivogen.com/pvitro2-luciaseap>

For research use only

Version 20H17-MM

PRODUCT INFORMATION

Contents:

- 20 µg of pVITRO2-blasti-Lucia/SEAP provided as lyophilized DNA
- 2 x 1 ml blasticidin at 10 mg/ml

Storage and stability:

- Product is shipped at room temperature.
- Upon receipt, store lyophilized DNA at -20°C.
- Resuspended DNA should be stored at -20°C.
- Store blasticidin at 4°C or -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

pVITRO is a family of vectors with improved features. pVITRO plasmids allow the co-expression of two or more genes from two different transcription units. pVITRO plasmids can be stably transfected in mammalian cells and are expressed at high levels.

pVITRO2-Lucia/SEAP contains the Lucia luciferase and SEAP reporter genes. pVITRO2-Lucia/SEAP can be used as a control vector.

pVITRO2-Lucia/SEAP also can be used for cloning of open reading frames (ORF). Both reporter genes are flanked by unique sites (BspH I/Avr II for Lucia luciferase and Nco I/Nhe I for SEAP) that allow for convenient cloning of ORFs.

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 water. Store resuspended plasmid at -20°C.

Plasmid amplification and cloning:

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

Blasticidin usage

Blasticidin should be used at 25-100 µg/ml in bacteria and 1-30 µg/ml in mammalian cells. Blasticidin is supplied at 10 mg/ml in HEPES buffer.

PLASMID FEATURES

- **hFerH and hFerL composite promoters:** Ferritin is a 24 subunit protein composed of two subunit types, termed H (heavy) and L (light), which perform complementary functions in the protein. Ferritin is ubiquitously expressed. Its synthesis is highly regulated by the iron status of the cell. The iron regulation is achieved at the translational level through the interaction between the iron-responsive element (IRE), located in the 5' untranslated region (5'UTR) of the ferritin mRNAs, and the iron regulatory protein¹. To eliminate the iron regulation of the ferritin promoters, the 5'UTR of FerH and FerL have been replaced by the 5'UTR of the mouse and chimpanzee elongation factor 1 (EF1) genes, respectively.

- **SV40 enhancer** which is comprised of a 72-base-pair repeat allows the enhancement of gene expression in a large host range. The enhancement varies from 2-fold in non-permissive cells to 20-fold in permissive cells. Furthermore, the SV40 enhancer is able to direct nuclear localization of plasmids².

- **CMV enhancer:** The major immediate early enhancer of the human cytomegalovirus (HCMV) is composed of unique and repeated sequence motifs. The HCMV enhancer can substitute for the 72-bp repeats of SV40 and is several-fold more active than the SV40 enhancer³.

- **pMB1 Ori** is a minimal *E. coli* origin of replication to limit vector size, but with the same activity as the longer Ori.

- **Lucia luciferase** is a synthetic CpG-free gene that codes for a secreted coelenterazine-utilizing luciferase. ORF size (from the ATG to the stop codon): 634 bp. Lucia luciferase activity can be evaluated using QUANTI-Luc™, an assay reagent containing all the components required to quantitatively measure the activity of Lucia luciferase and other coelenterazine-utilizing luciferases.

- **FMDV IRES:** The internal ribosome entry site of the Foot and Mouth Disease Virus enables the translation of two open reading frames from one mRNA with high levels of expression⁴.

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

- **Blasti:** Resistance to blasticidin is conferred by the *bsr* gene from *Bacillus cereus*. In bacteria, *bsr* is expressed from the constitutive *E. coli* EM7 promoter. In mammalian cells, *bsr* is transcribed from the hFerH/mEF1α promoter as a polycistronic mRNA and translated via the FMDV IRES.

- **EF1 pAn** is a strong polyadenylation signal. InvivoGen uses a sequence starting after the stop codon of the EF1 cDNA and finishing after a bent structure rich in GT.

- **SEAP** is a secreted form of human embryonic alkaline phosphatase. Unlike endogenous alkaline phosphatases, SEAP is extremely heat stable and resistant to the inhibitor L-homoarginine. It catalyses the hydrolysis of pNitrophenyl phosphate (pNpp) producing a yellow end product. SEAP expression can be readily quantified by collecting samples of culture medium and measuring the hydrolysis of pNpp with a spectrophotometer at 405 nm. SEAP activity that can be readily assessed qualitatively and quantitatively using HEK-Blue™ Detection or QUANTI-Blue™.

- **SV40 pAn:** The Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA. The efficiency of this signal was first described by Carswell *et al.*⁵

1. Kim DW. *et al.*, 1990. Use of the human elongation factor 1α promoter as a versatile and efficient expression system Gene 91:217-23. 2. Dean DA. *et al.*, 1999. Sequence requirements for plasmid nuclear import. Exp. Cell. Res. 253:713-22. 3. Boshart M. *et al.*, 1985. A very strong enhancer is located upstream of an immediate early gene of human cytomegalovirus. Cell 41:521-30. 4. Ramesh N. *et al.*, 1996. High-titer bicistronic retroviral vectors employing foot-and-mouth disease virus internal ribosome entry site. Nucleic Acids Res. 24:2697-700. 5. Carswell S. & Alwine JC. ,1989. Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. Mol. Cell Biol. 9: 4248-58.

TECHNICAL SUPPORT

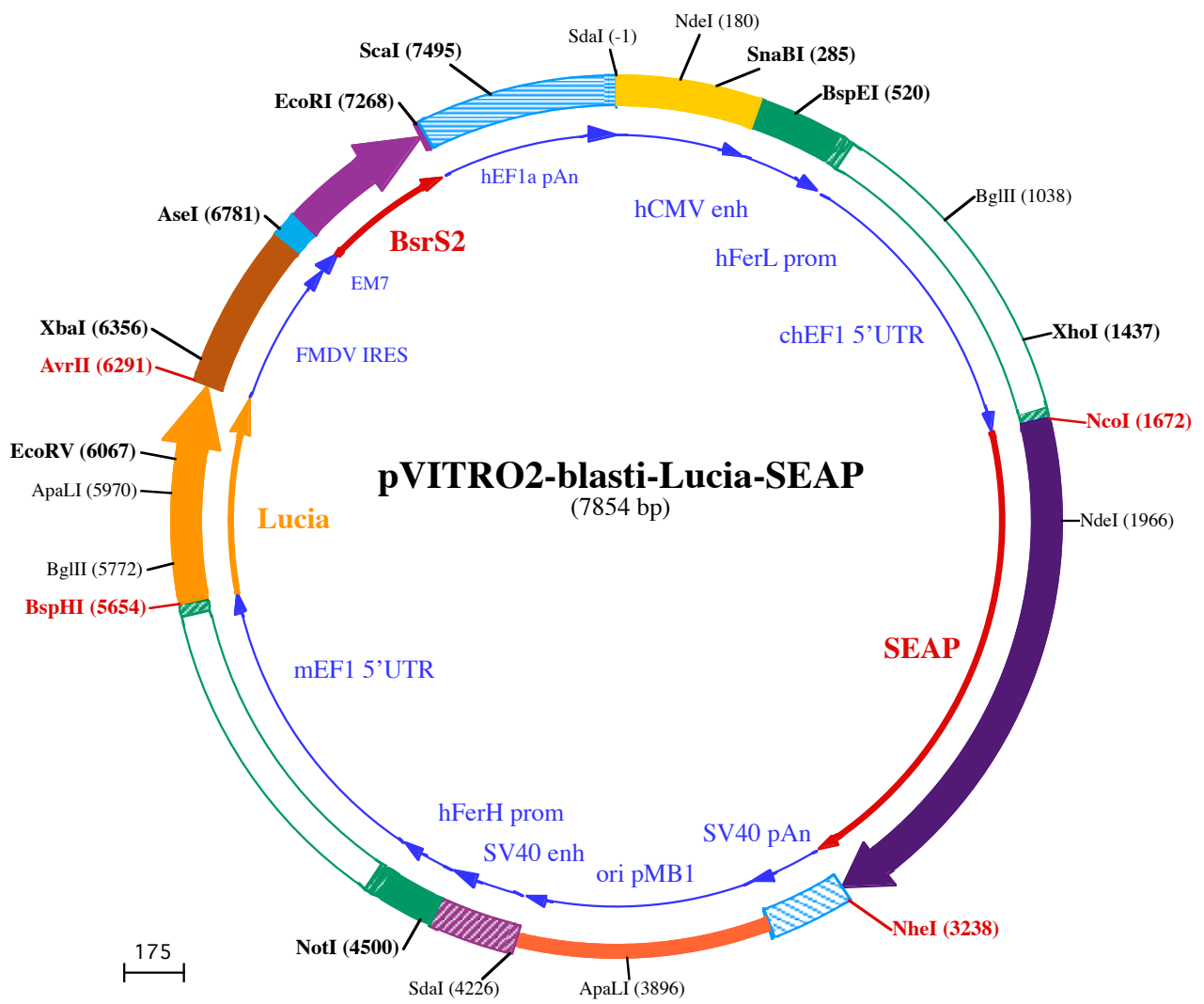
InvivoGen USA (Toll-Free): 888-457-5873

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



SdaI (-1)
1 CCTGCAGGCGTTACATAACTTACGGTAAATGGCCCGCTGGCTGACCGCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAA

NdeI (180)
101 CGCCAATAGGGACTTTCATTGACGTCAATGGGTGAGATTTACGGTAAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCC

SnaBI (285)
201 TATTGACGTCAATGACGGTAAATGGCCCGCTGGCATTATGCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATC
301 GCTATTACCATGATGATGCGGTTTTGGCAGTACATCAATGGGCGTGGATAGCGGTTTGACTIONCACGGGGATTCCAAGTCTCCACCCATTGACGTCAATG
401 GGAGTTTGTGTTGACTAGTCAGGGCCCAACCCCAAGCCCACTTTCACAACACGCTGGCGCTACAGGCGCTGACTTCCCCTTGCTTTGGGCGGG

BspEI (520)
501 GGGCTGAGACTCTATGTGCTCCGATTGGTCAGGCACGGCTTCGGCCCGCTCTGCCACCGAGATTGGCCGCTAGGCCTCCCGAGCGCCTGCC
601 TCCGAGGGCCGGCGACCATAAAGAAGCCCGCTAGCCACGTCCTCCGCGAGTTCGGCGGTCCCGGGTCTGTCTCAAGCTTGGCCGAGAACACAGg
701 taagtgcggtgtgtgttccccgggctgaccttttacgggttatggccttgcgtgccttgaattacttccatgcccctggctgcagtagctgattc
801 ttgatcccgagcttcgggttgaagtgggtgggagagttcaggccttgcgttaaggagcccttcgctcgtgcttgagttgagcctggcttggcg
901 ctggggccgctgctaactggtggcaccttcgctcgtcgtcgtcttctgctaagtctctagccatttaaaatgttataaccagctgcgag

BglIII (1038)
1001 cttttttctggcgagatagcttctaataatcgggccaagatctgcacactggtatcttgggttttggggccgcccggcgagccgggcccgtgcgtccc
1101 agcgacatgttcggcgaggccgggctgagcgcggccaccgagaatcggaggggtagctctaaactggccgctgctctggtgcctggcctcgc
1201 gccccgctgtatgccccccctgggggcaaggctggccggtggcaccagttgctgagcggaaagatggccgcttcccggcctgctcagggagc
1301 tcaaaatggaggacgcccggcgccgggagagcgggctgagtcacccacacaaaggaaggccttctctctcatccgctcgtctcatgtgactcca

XhoI (1437)
1401 cggagtaccggcgccgtccaggcacctcgattagttctcagcttttggagtacgctcgtcttaggttggggggaggggtttatgcatggagtttcc
1501 ccacactgagtggtggagactgaagagttaggccagcttggcacttgatgtaattctccttggaaattgcccctttttaggttggatcttgcctcattc

NcoI (1672)
1601 tcaagcctcagacagtggttcaagttttttcttccatttcagGTGTCGTGAAAACCTACCCCTAAAAGCCACCATGGTCTGGGGCCCTGCATGCTGT

1 M V L G P C M L L
1701 GCTGCTGCTGCTGGCCCTGAGGCTACAGCTCTCCCTGGGCATCATCCAGTTGAGGAGGAGAACCCTGGACTTCTGGAACCGCAGGCGAGCCGAGGCC
9 L L L L L G L R L Q L S L G I I P V E E E N P D F W N R E A A E A
1801 CTGGTGCCCAAGAAGCTGCAGCTGCACAGACAGCCCAAGAACCTCATCATCTTCTGGGCGATGGGATGGGGGTGTCTACGGTGACAGTGCACA
43 L G A A K K L Q P A Q T A A K N L I I F L G D G M G V S T V T A A

NdeI (1966)
1901 GGATCTAAAAGGGCAGAAGAAGGACAACTGGGGCTGAGATACCCTGGCTATGGACCGCTTCCCATATGTGGCTCTGTCCAAGACATAATGTAGA
76 R I L K G Q K K D K L G P E I P L A M D R F P Y V A L S K T Y N V D
2001 CAAACATGTGCCAGACAGTGGAGCCACAGCCACGGCCTACTGTGGGGGTCAAGGGCAACTTCCAGACCATGGCTTGTGAGTGCAGCCGCCCGCTTTAAC
109 K H V P D S G A T A T A Y L C G V K G N F Q T I G L S A A A R F N
2101 CAGTGAACACGACAGCGGGCAACGAGGTCTCTCCGTGATGAATCGGGCAAGAAAGCAGGGAAGTCAAGTGGGAGTGGTAACCCACACGAGTGCAGC
143 Q C N T T R G N E V I S V M N R A K K A G K S V G V V T T T R V Q
2201 ACGCTCGCCAGCCGACCTACGCCACACGGTGAACCGCACTGGTACTCGGACCGCCGACGTGCCTGCCTCGGCCCGCAGGAGGGTGCAGGACAT
176 H A S P A G T Y A H T V N R N W Y S D A D V P A S A R Q E G C Q D I
2301 CGTACGAGCTCATCTCAACATGGACATTGATGTGATCTGGTGGAGGCCGAAAGTACATGTTTCGCATGGGAACCCAGACCCTGAGTACCCAGAT
209 A T Q L I S N M D I D V I L G G R K Y M F R M G T P D P E Y P D
2401 GACTACGCCAAGTGGGACCAGGCTGAGCGGAAGAATCTGGTGCAGGAATGGCTGGCGAAGCGCCAGGGTCCCGGTATGTGTGAACCCGACTGAGC
243 D Y S Q G G T R L D G K N L V Q E W L A K R Q G A R Y V W N R T E
2501 TCATGCAGGCTTCCCTGGACCCGTCTGTGACCCATCTCATGGGTCTCTTTGAGCCTGGAGACATGAAATACGAGATCCACCGAGACTCCACACTGGACCC
276 L M Q A S L D P S V T H L M G L F E P G D M K Y E I H R D S T L D P
2601 CTCCTGATGGAGATGACAGAGGCTGCCCTGCGCTGCTGAGCAGGAACCCCGGGCTTCTTCTCTTCTGTTGGAGGGTGGTGCATCGACCACGGTCA
309 S L M E M T E A A L R L L S R N P R G F F L F V E G G R I D H G H
2701 CACGAAAGCAGGGCTTACCGGCACTGACTGAGACGATCATGTTTCGACGACGCCATTGAGAGGGCGGGCCAGCTCACGAGGAGGACACGCTGAGCC
343 H E S R A Y R A L T E T I M F D D A I E R A G Q L T S E E D T L S
2801 TCGTCACTGCCGACCACTCCACGCTTCTCTTCCGGAGGCTACCCCTCGGAGGGAGCTCCATCTTGGGCTGGCCCTGGCAAGCCCGGACAGGAA
376 L V T A D H S H V F S F G G Y P L R G S S I F G L A P G K A R D R K

2901 GGCCTACACGGTCTCTATACGAAACGGTCCAGGCTATGTGCTCAAGGACGGCGCCCGCGGATGTTACCAGAGCGAGAGCGGGAGCCCGAGTAT

409▶ A Y T V L L Y G N G P G Y V L K D G A R P D V T E S E S G S P E Y

3001 CGGCAGCAGTCAGCAGTGCCTTGACGAAGAGACCCACGAGGCGAGGACGTGGCGGTGTTCCGCGCGGCCGAGGCGCACCTGGTTCACGGCGTGC

443▶ R Q Q S A V P L D E E T H A G E D V A V F A R G P Q A H L V H G V

3101 AGGAGCAGACCTTCATAGCGCACGTCATGGCTTCGCGCGCTGCCTGAGGCTACACCGCTGCGACCTGGCGCCCGCCGGCACCACCGACCGCCG

476▶ Q E Q T F I A H V M A F A A C L E P Y T A C D L A P P A G T T D A A

NheI (3238)

3201 GCACCCGGGGCGTCCCGTCCAAGCGTCTGGATTGAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCA

509▶ H P G R S R S K R L D •

3301 GTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAGTTAAACAACAACATTCATTCAT

3401 TTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAAACCTCTACAATGTGGTATGGAATGTTAATTAAGTACCATGACCAA

3501 AATCCCTTAACGTGAGTTTTTCGTTCCACTGAGCGTCAGACCCGTAGAAAAGATCAAAGGATCTTCTTGAGATCCTTTTTTTCGCGGTAATCTGCTGC

3601 TTGCAACAAAAAACCCCGCTACCAGCGGTGGTTGTTTGGCGGATCAAGAGCTACCAACTCTTTTTCCGAAGGTAAGTGGCTTCAGCAGAGCGCAGA

3701 TACCAAATACTGTTCTTCTAGTGTAGCCGTAGTTAGGCCACCACTCAAGAACTCTGTAGCACCGCTACATACCTCGCTCTGCTAATCCTGTTACCAGT

ApaI (3896)

3801 GGCTGCTGCCAGTGGCGATAAGTCGTGCTTACCGGTTGGACTCAAGACGATAGTTACCGGATAAGGCGCAGCGGTGGGCTGAACGGGGGTTCTGTGC

3901 ACACAGCCAGCTTGGAGCGAACGACCTACCCGAACTGAGATACCTACAGCGTGTGATGAGAAAGCGCCAGCTTCCGAAGGGAGAAAGCGGACA

4001 GGTATCCGGTAAGCGCGAGGTCGGAACAGGAGAGCGCACGAGGAGCTTCCAGGGGAAACGCCTGGTATCTTTATAGTCTGTCGGGTTTCGCACCT

4101 CTGACTTGAGCGTCGATTTTTGTGATGCTGTCAGGGGGCGGAGCCTATGAAAAACGCCAGCAACGCGGCTTTTTACGGTTCCTGGCCTTTTGTGG

SdaI (4226)

4201 CCTTTTGCTCACATGTTCTTAATTAACCTGCAGGGCTGAAATAACCTCTGAAAGAGGAACCTGGTTAGGTACCTTCTGAGGCTGAAAGAACCAGCTGTG

4301 GAATGTGTGTCAGTTAGGGTGTGAAAGTCCCGAGGCTCCCGAGCAGGAGGATGCAAAAGCATGCATCTCAATTAGTCAGCAACCAGGTGTGAAAG

4401 TCCCCAGGCTCCCGAGCAGGAGGATGCAAAAGCATGCATCTCAATTAGTCAGCAACCATAGTCCCACTAGTTCCGCCAGAGCGCGAGGGCTCCA

NotI (4500)

4501 GCGGCCGCCCTCCCGCAGCAGGGGCGGGTCCCGCGCCACCGAAGGAGCGGGCTCGGGCGGGCGGCTGATTGGCGGGGCGGGCTGACGCC

4601 GACGCGGTATAAGAGACCAAGCGACCCGAGGGCAGAGCTTCTTCGCCAAGCTTGGCGTCAGAACGAGGTGAGGGGCGGGTGTGGTTCGCGGG

4701 GCCGCGAGCTGGAGTCTGCTCCGAGCGGGCGGGCCCGCTGCTGCTCGCGGGGATTAGCTGCGAGCATTCCCGCTTCGAGTTGCGGGCGCGCGGG

4801 AGGCAGAGTGCAGGCTAGCGCAACCCGTAGCCTCGCTCGTGTCCGGCTTGGAGCTAGCGTGGTGTCCGCGCCCGCCGCGTGTACTCCGGCC

4901 GCACTCTGGCTTTTTTTTTTTTGTGTTGTTGCCCTGCTGCTTCGATTGCGGTTGAGCAATAGGGGTAACAAGGGAGGGTGCGGGCTTGTGCGCC

5001 CGGAGCCCGAGAGGTGATGTTGGGAGGAATGGAGGGACAGGAGTGGCGCTGGGGCCCGCCGCTTCGGAGCACATGTCGACGCCACCTGGATGG

5101 GCGGAGGCTGGGTTTTTCCGAAGCAACAGGCTGGGTTAGCGTGCCGAGGCCATGTGGCCCGCAGCACCCGGCAGATCTGGCTTGGCGGCGCCGG

5201 TTGCCCTGCTCCCTAAGAGGTGAGCCATCCCGTCCGCGCAGTGGCTGCTGGAAGATGGCCGCTCCCGGCCCTGTTGCAAGGAGCTCAAAA

5301 TGGAGGACGCGGAGCCCGGTGGAGCGGGGGTGGTACCCACACAAAGGAGGGCTGGTCCCTACCGGCTGCTGCTTCTGTGACCCGTGGT

5401 CCTATCGGCCAATAGTCACCTCGGGCTTTGAGCAGGCTAGTCGCGGGGGGAGGGATGTAATGGCGTTGGAGTTTGTTCACATTTGGTGGGTG

5501 GAGACTAGTCAGGCCAGCCTGGCGTGAAGTCATTTTTGGAATTTGTCCTTGGATTTTGGAGCGGAGCTAATTCGGGCTTCTTAGCGGTTCAAAGG

BspHI (5654)

5601 TATCTTTAAACCTTTTTTGTGTTGTGAAAACCACCGTAATCAAAGCAATCATGATGGAATCAAGGTGCTGTTGCCCTCATCTGATTGCTGT

1▶ M M E I K V L F A L I C I A V

BglIII (5772)

5701 TGCTGAGGCAAAACCCACTGAAATCAATGAAGACCTCAATATAGTGTGCTGGCCCTCAACTTTGCCACCACAGATCTTGAGACTGACCTGTTCAACCAAC

15▶ 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 N

5801 TGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGCAAGTGCCTGGCAAAAACCTCCCGCAGATGCTCTGA

49▶ 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 L

ApaI (5970)

5901 GGGAGCTGGAGCCAATGCCAGAAGGGCTGGTGCACAAGAGGCTGCTCATTGCTCTCCACATTAAGTGACCCCTAAGATGAAGAAATTTATCCC

82▶ 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 P

EcoRV (6067)

6001 TGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCTGGCTCAAGGATAAG

115▶ 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 K
6101 GAGCCACTGGACCAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGACCACTGGCTGTCTGAAGGGCCTTGCCAATGTCCAGTGTCTGACCTCTGA

149▶ 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 L

6201 AGAAGTGGCTTCCCCAGAGGTGTACCACCTTTTCCAGCAAGATTACAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGATGATACCTAGGAGC **AvrII (6291)**

182▶ 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 •

XbaI (6356)

6301 AGGTTTCCCAATGACACAAAACGTGCAACTTGAAGTCCGCTGGTCTTTCCAGGTCTAGAGGGTAACACTTTGTAAGTGGCTTGGCTCCACGCTCGA

6401 TCCACTGGCGAGTGTAGTAACAGCACTGTTGCTTGTAGCGGAGCATGACGGCCGTGGAACTCCTCCTTGGTAACAAGGACCCACGGGGCCAAAAGCC

6501 ACGCCACACGGGCCGTGATGTGTCAACCCAGCACGGCGACTTACTGCGAAACCACTTTAAAGTGACATTGAAACTGGTACCCACACTGTTGA

6601 CAGGCTAAGGATGCCCTTCAAGTACCCGAGGTAAACACGCGACACTCGGGATCTGAGAAGGGACTGGGGCTTCTATAAAAGCGCTCGGTTTAAAAAGCT

AseI (6781)

6701 TCTATGCCTGAATAGGTGACCGGAGGTCGGCACCTTTCCTTTGCAATTACTGACCTATGAATACAA **CTGACTGTTTGACAATTAATCATCGGCATAGTA**

6801 **TATCGGCATAGTATAATACGACTCACTATAGGAGGGCCACCATGAAGACCTTCAACATCTCTCAGCAGGATCTGGAGCTGGTGGAGGTCGCCACTGAGAA**

1▶ M K T F N I S Q Q D L E L V E V A T E K

6901 GATCACCATGCTCTATGAGGACAACAAGCACCATGTGGGGCGGCCATCAGGACCAAGACTGGGGAGATCATCTCTGCTGCCACATTGAGGCCTACATT

20▶ I T M L Y E D N K H H V G A A I R T K T G E I I S A V H I E A Y I
7001 GGCAGGGTCACTGTCTGTCTGAAGCATTGCCATTGGTCTGCTGTGAGCAACGGGCAGAAGGACTTTGACACCATTGTGGCTGCAGGCCACCCCTACT

54▶ G R V T V C A E A I A I G S A V S N G Q K D F D T I V A V R H P Y
7101 CTGATGAGGTGGACAGATCCATCAGGGTGGTCAGCCCTGTGGCATGTGACAGAGCTCATCTCTGACTATGCTCCTGACTGCTTTGTGCTCATTGAGAT

87▶ S D E V D R S I R V V S P C G M C R E L I S D Y A P D C F V L I E M

EcoRI (7268)

7201 GAATGGCAAGCTGGTCAAAACACCATTGAGGAACATCCCCCTCAAGTACACCAGGAACATAACCTGAATTGCTAGG **ATTATCCCTAATACCTGCCA**

120▶ N G K L V K T T I E E L I P L K Y T R N •
7301 CCCCCTCTTAATCAGTGGTGAAGAACGGTCTCAGAACTGTTGTTCAATTGGCCATTTAAGTTTAGTAGTAAAAGACTGGTTAATGATAACAATGCA

ScaI (7495)

7401 TCGTAAAACCTTCAGAAGGAAAGGAGAATGTTTTGTGACCACCTTGGTTTTCTTTTTGCGTGTGGCAGTTTTAAGTTATTAGTTTTAAAATCAGTAC

7501 TTTTTAATGAAAACAATTGACCAAAAATTTGTACAGAATTTGAGACCCATTAATAAGTAAATGAGAAACCTGTGTTCCTTTGGTCAACACCGA

7601 GACATTTAGGTGAAAGACATCTAATTCTGTTTTACGAATCTGGAACTTCTTAAAAATGTAATCTTGAGTTAACACTTCTGGGTGGAGAATAGGGTTG

7701 TTTTCCCCCACAATTTGGAAGGGGAAGGAATATCATTTAAAGCTATGGGAGGGTGTCTTTGATTACAACACTGGAGAGAAATGCAGCATGTTGCTGAT

7801 TGCCTGCACTAAAACAGGCCAAAAACTGAGTCTTGGGTGCATAGAAAAGCTG