

STOP

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

Important Limited Use License information for pVITRO1-neo-Lucia/SEAP

The purchase of the pVITRO1-neo-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 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



pVITRO1-neo-Lucia/SEAP

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

Catalog code: pvitro1-nlucsp

<https://www.invivogen.com/pvitro1-luciaseap>

For research use only

Version 19L17-MM

PRODUCT INFORMATION

Contents:

- 20 µg of pVITRO1-neo-Lucia/SEAP provided as lyophilized DNA

Storage and stability:

- Product is shipped at room temperature.
- Lyophilized DNA is stable for 3 months at -20°C.
- Resuspended DNA is stable more than one year at -20°C.

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 new 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.

pVITRO1-Lucia/SEAP contains the reporter genes Lucia luciferase and SEAP and can be used as a control vector.

pVITRO1-Lucia/SEAP also can be used for cloning of open reading frames (ORF). Both reporter genes are flanked by unique sites (*NcoI*/*AvrII* for Lucia luciferase and *BspHI*/*NheI* for SEAP) that allow for convenient cloning of ORF's

PLASMID FEATURES

- **rEF1 and mEF1 prom:** pVITRO1-neo-GFP/LacZ plasmid carries two elongation factor 1 alpha (EF-1α) promoters, from rat and mouse origins. Similarly to their human counterpart¹, both promoters display a strong activity that yield similar levels of expression. EF-1α promoters are expressed at high levels in all cell cycles and lower levels during G0 phase. EF-1α promoters are also non-tissue specific; they are highly expressed in all cell types.
- **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.
- **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.*⁴
- **pMB1 Ori** is a minimal *E. coli* origin of replication to limit vector size, but with the same activity as the longer Ori.
- **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 expression⁵.
- **EM7** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*.

- **Neo:** The neo gene from Tn5 encodes an aminoglycoside 3'-phosphotransferase (3' APH II) that confers resistance to the antibiotics kanamycin in bacteria and G418 in mammalian cells. In bacteria, *neo* is expressed from the constitutive *E. coli* EM7 promoter. In mammalian cells, *neo* is transcribed from the rat EF-1α 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.

- **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.

- **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™.

1. Kim DW. *et al.*, 1990. Use of the human elongation factor 1α promoter as a versatile and efficient expression system Gene 91(2):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(2):521-30. 4. 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. 5. Ramesh N. *et al.*, 1996. High-iter bicistronic retroviral vectors employing foot-and-mouth disease virus internal ribosome entry site. Nucleic Acids Res. 24(14):2697-700.

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 other commonly used laboratory *E. coli* strains, such as DH5α.

Bacterial antibiotic selection

Kanamycin (not provided) is normally used for *E. coli* at a final concentration of 50 µg/ml in liquid or solid media.

Mammalian antibiotic selection

G418 is normally used at a concentration of 400 µg/ml. However, the optimal concentration needs to be determined for your cells.

RELATED PRODUCTS

Product	Description	Cat. Code
ChemiComp GT116 cells	Competent <i>E. coli</i> cells	gt116-11
G418	Selection antibiotic	ant-gn-1

TECHNICAL SUPPORT

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

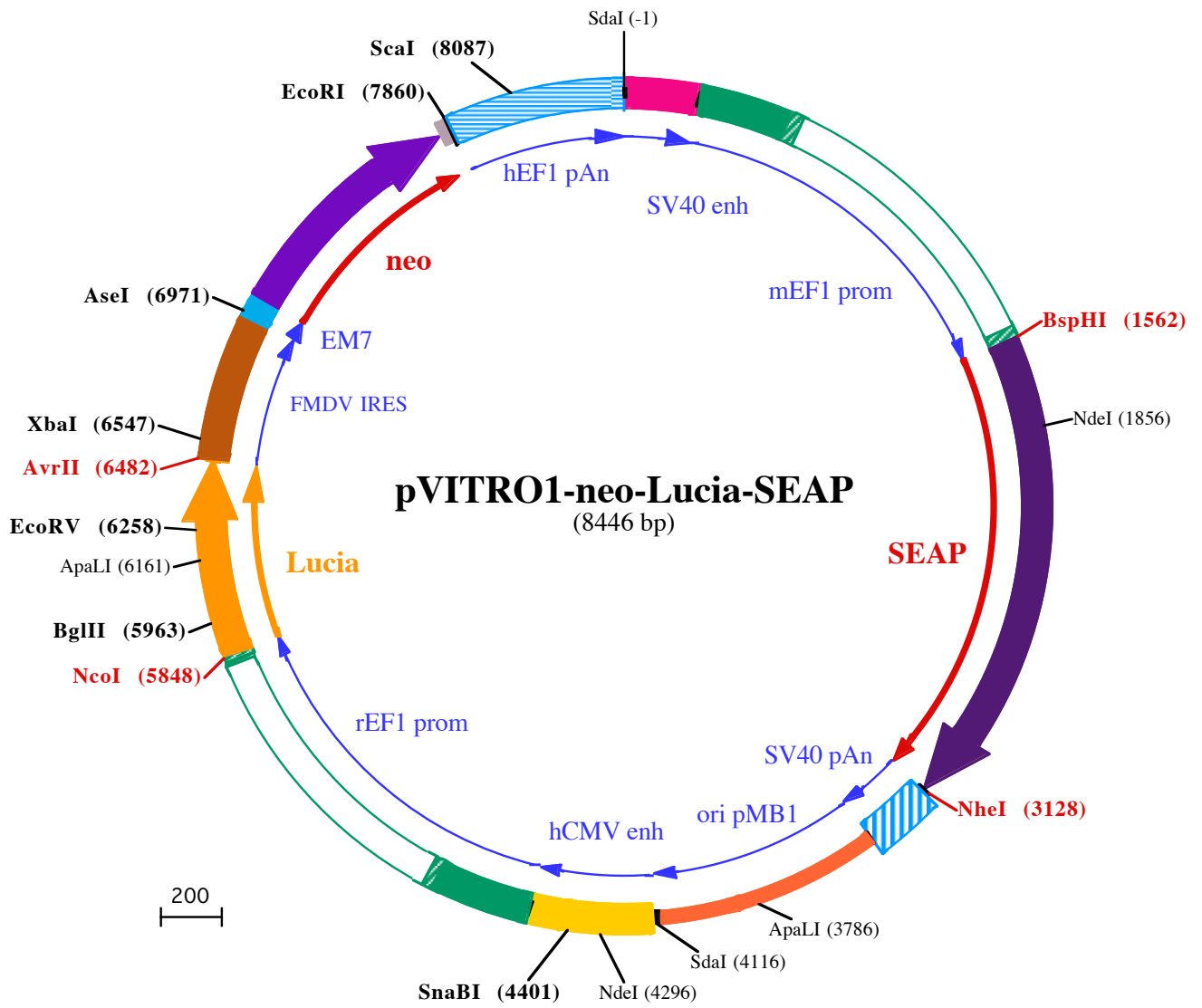
InvivoGen USA (International): +1 (858) 457-5873

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

InvivoGen Hong Kong: +852 3622-3480

E-mail: info@invivogen.com

 **InvivoGen**
www.invivogen.com



SdaI (-1)
1 CCTGCAGGGCCTGAAATAACCTCTGAAAGAGGAACCTTGGTTAGGTACCTTCTGAGGCGAAAGAACCAGCTGTGAATGTGTTCAGTTAGGGTGTGGAA

101 AGTCCCCAGGCTCCCCAGCAGGCAGAAGTATGCAAAGCATGCATCTCAATTAGTCAGCAACCAGGTGTGAAAGTCCCCAGGCTCCCCAGCAGGCAGAAG

201 TATGCAAAGCATGCATCTCAATTAGTCAGCAACCATAGTCCACTAGTGGAGCCGAGAGTAATTCATACAAAAGGAGGGATCGCTTCGCAAGGGGAGAG

301 CCCAGGGACCGTCCCTAAATTTCTACAGACCCAAATCCCTGTAGCCGCCACGACAGCGGAGGAGCATGCGCTCAGGGCTGAGCGCGGGGAGAGCAGA

401 GCACACAAGCTCATAGACCCTGGTCGTGGGGGGAGGACCGGGGAGCTGGCGGGGCAAACCTGGGAAAGCGGTGTCGTGTCTGCTCCGCCCTTTCC

501 CGAGGGTGGGGGAGAACGGTATATAAGTGGCGCAGTCGCTTGGACGTTCTTTTTTCGCAACGGGTTTGCCGTCAGAACCGAGGTGAGGGGCGGGTGTGGC

601 TTCCGCGGGCCCGCAGCTGGAGGTCCTGCTCCGAGCGGGCCGGCCCGCTGTCGTCGCGGGGATTAGCTGCGAGCATTCCGCTTCGAGTTGCGGGC

701 GCGCGGGAGGCAGAGTGCAGGCGCTAGCGGCAACCCGTAGCCTCGCTCGTGTCCGGCTTGGAGCCTAGCGTGGTGTCCGCGCCGCCCGCGCTGCTA

801 CTCCGGCCGACTCTGCTTTTTTTTTTTTTTTTGTGTTGTTGCCCTGCTGCCTTCGATTGCCGTTACGAATAGGGGCTAACAAAGGGAGGGTGGCGGGCT

901 TGCTCGCCCGAGCCCGAGAGGTGATGTTGGGGAGGAATGGAGGGACAGAGTGGCGGCTGGGGCCCGCCGCTTCGAGACATGTCCGAGCCAC

1001 CTGGATGGGGCAGGCTGGGGTTTTTCCGAAGCAACCAGGCTGGGGTTAGCGTGCCGAGGCCATGTGGCCACGACCCGGCAGCATGTGGCTTGGCG

1101 GCGCCGCTTGCCTGCCTCCCTAACTAGGGTGAGGCCATCCGTCGGCCACCGTTGCGTGCCTGAAAGATGGCGCTCCCGGGCCCTGTTGCAAGGA

1201 GCTCAAATGGAGACCGCGCAGCCGGTGGAGCGGGGGTGGAGTACCCACACAAAGGAAGGGCCTGGTCCCTACCGGCTGCTGCTTCTGTGAC

1301 CCCGTGCTCTATCGCCGCAATAGTCACTCGGGCTTTTGGAGCAGGCTAGTCGCGCGGGGGAGGGATGTAATGGCGTTGGAGTTTGTTCACATTT

1401 GGTGGTGGAGACTAGTCAGGCCAGCCTGGCGCTGGAAGTATTTTTGGAATTTGTCCCTTGGATTTTGGCGGAGCTAATTTCTCGGGCTTCTTAGCGG

1501 TTCAAAGGTATCTTTTAAACCTTTTTTAGGTGTTGTGAAAACACCGCTAATTCAAAGCAATCATGATTCTGGGCGCTGCATGCTGCTGCTGCTGCTG

BspHI (1562)

1▶ M I L G P C M L L L L L L

1601 CTGCTGGGCGTGGGCTACAGCTCTCCCTGGGCATCATCCAGTTGAGGAGGAGAACCAGGACTTCTGGAACCGCAGGCGAGCCGAGGCGCTGGGTGCCG

13▶ 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 L G A

1701 CCAAGAAGCTGCAGCCTGCACAGACAGCCGCAAGAACCTCATCATCTTCTGGCGATGGGATGGGGGTGTCTACGGTGACAGCTGCCAGGATCCTAAA

46▶ 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 R I L K

NdeI (1856)

1801 AGGGCAGAAGAAGGACAAAACCTGGGCGCTGAGATACCCCTGGCTATGGACCGCTTCCCATATGTGGCTCTGTCCAAGACATAAATGTAGACAAAATGTG

79▶ 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 K H V

1901 CCAGACAGTGGAGCCACAGCCAGGCGCTACCTGTGGGGGTCAAGGGCAACTTCCAGACCATTTGGCTTGGTGCAGCGCCCGCTTAAACAGTGAACA

113▶ 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 Q C N

2001 CGACACGGCAACGAGGTGATCTCCGCTGATGAATCGGGCCAAGAAAGCAGGGAAGTCAAGTGGGAGTGGTAACCCACACGAGTGCAGCAGCGCTCGCC

146▶ 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 H A S P

2101 AGCCGGCACCTACGCCACAGGTAACCGCAACTGTTACTCGGACCGCAGCTGCCTGCCTCGGCCCGCAGGAGGGTGCAGGACATCGCTACGCGA

179▶ 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 A T Q

2201 CTCATCTCAACATGGACATTTGATGTGATCTGGGTGGAGGCCGAAAGTACATGTTTCGATGGGAACCCAGACCCCTGAGTACCAGATGACTACAGCC

213▶ L I S N M D I D V I L G G G R K Y M F R M G T P D P E Y P D D Y S

2301 AAGGTGGACAGGCTGGAGCGGAAGAACTGTTGTCAGGAATGGCTGGCGAAGCGCCAGGGTGCCTGGTATGTGTGAACCGCACGAGCTCATGACGGC

246▶ 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 L M Q A

2401 TTCCCTGGACCGTCTGTGACCATCTCATGGTCTCTTTGAGCTGGAGACATGAAATACGAGATCCACCAGACTCCACACTGGACCCCTCCCTGATG

279▶ 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 S L M

2501 GAGATGACAGAGGCTGCCCTGGCGCTGCTGAGCAGGAACCCCGGGCTTCTCTCTTCTGTTGGAGGGTGGTGCATCGACCAGGTCATCACGAAAGCA

313▶ 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 H E S

2601 GGGCTTACCGGCACTGACTGAGACGATCATGTTGACGACGCCATTGAGAGGGCGGGCAGCTCACCAGCGAGGAGACACGCTGAGCCTCGTCACTGC

346▶ 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 L V T A

2701 CGACCACTCCACGCTTCTCTTCCGAGGCTACCCCTGCGAGGGAGCTCCATCTTCCGGCTGGCCCTGGCAAGGCCGGGACAGGAAGGCTACAGC

379▶ 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 A Y T

2801 GTCCTCTATACGAAACGGTCCAGGCTATGTGCTCAAGGACGGCGCCCGGGATGTTACCGAGAGCGAGAGCGGGAGCCCGAGTATCGGACGAGT

413▶ 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 R Q Q

2901 CAGCAGTGGCCCTGGACGAAGAGACCCACGAGCGAGGAGCTGGCGGTGTTGCGCGCGGGCCCGAGGCGCACCTGGTTCACGGCGTGCAGGAGCAGAC

446▶ 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 Q E Q T

3001 CTTATAGCGCAGTGCATGGCTTCCGCGCTGCTGGAGCCCTACACCGCTGCGACCTGGCGCCCGCCGGCACCACCGACCGCGCCACCCGGG

479▶ 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 H P G

NheI (3128)

3101 CGTCCCGGTCGAAGCTGCTGGATTGAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAACCCACAACCTAGAATGCAGTGAAAAAAA

513▶ R S R S K R L D •

3201 TGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAACAAATTGCATTCAATTTATGTTTC

3301 AGGTTTCAGGGGAGGTTGGGGAGGTTTTTAAAGCAAGTAAACCTCTACAATATGGTATGAAATGTTAATTAACACTAGCCATGACCAAAATCCCTTAA

3401 CGTGAGTTTTCTGTTCCACTGAGCGTCAGACCCCGTAGAAAAGATCAAAGGATCTTCTTGGAGATCCTTTTTTCTGCGGTAATCTGCTGCTTGCACAAACA

3501 AAAAACCCACCGCTACCAGCGGTGGTTTGGTTTCCGGATCAAGAGCTACCAACTCTTTTCCGAAGTAACTGGCTTCAGCAGAGCGCAGATACCAATAC
3601 TGTTCTTCTAGTGTAGCCGTAGTTAGGCCACCACCTCAAGAACTCTGTAGCACCGCTACATACCTCGCTCTGTAATCTGTTACCAGTGGCTGTGCC
3701 AGTGGCGATAAGTCGTGTCTTACCAGGTTGGACTCAAGACGATAGTTACCAGATAAGGCGCAGCGGTCGGGCTGAACGGGGGGTTCTGTGCACACAGCCCA
3801 GCTTGGAGCGAACGACCTACACCGAACTGAGATACCTACAGCGTGAGCTATGAGAAAGCGCCACGCTTCCCAGAGGGAGAAAGCGGACAGGTATCCGGT
3901 AAGCGCGAGGGTCGGAACAGGAGAGCGCACGAGGGAGCTTCCAGGGGAAACGCTGGTATCTTTATAGTCTGTGGGTTTCGCCACCTCTGACTTGAG
4001 CGTCGATTTTTGTGATGCTCGTCAGGGGGCGGAGCCTATGAAAAACGCCAGCAACGCGGCTTTTTACGGTTCCTGGCCTTTTGTGGCCTTTTGTCTC
4101 ACATGTTCTTAATTAACCTGCAGGCGTTACATAACTTACGGTAAATGGCCCGCTGGCTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACG
4201 TATGTTCCCATAGTAACGCCAATAGGGACTTTCCATTGACGTCAATGGGTGGAGTATTTACGGTAACTGCCACTTGGCAGTACATCAAGTGTATCATA
4301 TGCCAAGTACGCCCTATTGACGTCAATGACGGTAAATGGCCCGCTGGCATTATGCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACAT
4401 CTACGTATTAGTCATCGCTATTACCATGATGATGCGGTTTTGGCAGTACATCAATGGGCGTGGATAGCGGTTTGACTCACGGGGATTTCCAAGTCTCCAC
4501 CCCATTGACGTCAATGGGAGTTTTGTTTTACTAGTGGAGCCGAGAGTAATTCATACAAAAGGAGGATCGCCTTCGAAGGGGAGAGCCAGGGACCGTC
4601 CCTAAATTCTCACAGACCCAAATCCCTGTAGCCGCCACGACAGCGCAGGAGCATGCGCCAGGGCTGAGCGCGGTAGATCAGAGCACACAAGCTCA
4701 CAGTCCCGCGGTTGGGGGAGGGGCGCTGAGCGGGGCGAGGAGCTGGCGCGGGCAAACCTGGAAAGTGGTGTCTGTGGTCCGCCCTCTT
4801 CCCGAGGGTGGGGGAGAACGGTATATAAGTCGCGTAGTCGCCTTGACGTTCTTTTTCGCAACGGGTTTGGCCTCAGAACGCAgtagtgagggtg
4901 gcttccgaggccccggagctggagccctgctctgagcgggccccgctgatatgagtgctgctcgcagggttagctgtgagcattcccacttcgagt
5001 ggcgggctgaggggggagagtgagagcctagcggaacccccctagcctcgcctcgtgctcggcttaggcctagcgtggtgctccgcccgcgctg
5101 cactccgcccactatgctgtttttgctctgctgctcctcattgctcctcagcagcatgggctaacaaggagggtggtgggctcactcttaaggag
5201 cccatgaagcttacggttgataggaatggaaggcaggaggggcactggggcccgccttcggagcacatgtccgacgccacctggatggggcgag
5301 gcctgtggctttccgaagcaatcgggctgagtttagcctacctgggcatgtggccctagcactgggcacggctggtcctggcgggtgcccgttcctt
5401 gcctccaacaagggtgagggcctccgcccggcaccagttgctgctgagcggaaagatggcgctccggggccctgttgcaaggagctcaaaatggagga
5501 cgcggcagcccgggtgagcgggctgagtcaccacacaaggaagaggccttgcctcgcggcggctgcttctgtgaccccgtggtctatcgg
5601 ccgcatagtcacctcgggcttctcttgagcaccgctcgtcgcggcgggggaggggatctaaggcgttgagttgttcacatttgggtgggagact
5701 agtcaggccagcctggcgtggaagctattcttgaatttcccccttgagttggagcgagctaattctcaagcctcttagcgttcaaaggatattt
5801 ctaaaccggtttccagGTGTTGTGAAAGCCACCGCTAATTCAAAGCAACcATGGAAATCAAGGTGCTGTTTCCCTCATCTGTATTGCTGTTGCTGAGGC
1▶ M E I K V L F A L I C I A V A E A
5901 AAAACCCACTGAAATCAATGAAGACCTCAATATAGCTGTGGCCTCCAACCTTGGCCACCACAGATCTTGAGACTGACCTGTTACCAACTGGGAGACC
17▶ 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 W E T
6001 ATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACTCCCCCAGATGCTCTGAGGGAGCTGG
51▶ 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 R E L
6101 AGGCCAATGCCAGAAGGCTGGTTGCACAAGAGGCTGCCTCATTTGCCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTG
84▶ 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 G R C
6201 CCACACTTATGAAGGTGAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCTGGCTTCAAGGATAAGGAGCCACTG
117▶ 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 E P L
6301 GACCAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGACCACTGGCTGTCTGAAAGGCTTGCATGTCAGTGTCTGACCTCCTGAAGAAGTGGC
151▶ 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 K K W
6401 TTCCCCAGAGGTGACCACTTTTCCAGCAAGATTAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGATGATACCTAGGAGCAGGTTTCCC
184▶ 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 •
6501 CAATGACACAAAACGTGCAACTTGAAGCTCCGCTGGTCTTCCAGGTCTAGAGGGGTAACACTTTGACTGCGTTTGGCTCCAGCTCGATCCACTGGC
6601 GAGTGTAGTAACAGCACTGTTGCTTCTGAGCGGAGCATGACGGCCGTGGGAACTCCTCCTTGGTAACAAGGACCCAGGGGCCAAAAGCCAGCCCA
6701 CGGGCCCGTCACTGTGTCAACCCAGCAGCGGACTTTACTGCGAAACCACTTTAAAGTGCATTGAAAAGTGGTACCCACACACTGGTACAGGCTAAG
6801 GATGCCCTTCAAGTACCCGAGGTAACACGCGACACTCGGGATCTGAGAAGGGGACTGGGCTTCTATAAAGCGCTCGGTTAAAAAGCTTCTATGCCT
6901 GAATAGGTGACCGGAGGTGGCACCTTCTTTGCAATTACTGACCTATGAATACACTGACTGTTGACAATTAATCATCGGCATAGTATATCGGCAT
7000 AGTATAATACGACTCACTATAGGAGGGCCACCATGATTGAACAAGATGAGTTGACGCGAGGTTCTCCGGCCGCTGGGTGGAGAGGCTATTGCGCTATGA
1▶ M I E Q D G L H A G S P A A W V E R L F G Y D

7100 CTGGGCACAACAGACAATCGGCTGCTCTGATGCCCGGTGTTCCGGCTGTCAGCGCAGGGCGCCCGTTCTTTTGTCAAGACCGACCTGTCCGGTGCC
23▶ W A Q Q T I G C S D A A V F R L S A Q G R P V L F V K T D L S G A
7200 CTGAATGAAGTGAAGACGAGCGAGCGCGGCTATCGTGGCTGGCCACGACGGCGTTCCTTGCGCAGCTGTGCTCGACGTTGCTACTGAAGCGGAAGG
57▶ L N E L Q D E A A R L S W L A T T G V P C A A V L D V V T E A G R
7300 ACTGGCTGCTATTGGCGAAGTGGCGGGCAGGATCTCTGTCATCTCACCTTGTCTCTGCCGAGAAAGTATCCATCATGGCTGATGCAATGCGGCGGCT
90▶ D W L L L G E V P G Q D L L S S H L A P A E K V S I M A D A M R R L
7400 GCATACGCTTGATCCGGCTACTGCCATTGACCACCAAGCGAAACATCGCATCGAGCGAGCACGTACTCGGATGGAAGCCGGTCTTGTGATCAGGAT
123▶ H T L D P A T C P F D H Q A K H R I E R A R T R M E A G L V D Q D
7500 GATCTGGACGAAGAGCATCAGGGCTCGCGCCAGCCGAAGTTCGCCAGGCTCAAGGCGAGCATGCCCGAGCGGAGGATCTCGTGTGACACATGGCG
157▶ D L D E E H Q G L A P A E L F A R L K A S M P D G E D L V V T H G
7600 ATGCTGCTTGCCGAATATCATGGTGGAAAATGGCCGCTTTCTGGATTCATCGACTGTGGCCGGCTGGGTGTGGCGGACCCGCTATCAGGACATAGCGTT
190▶ D A C L P N I M V E N G R F S G F I D C G R L G V A D R Y Q D I A L
7700 GGCTACCGTGATATTGCTGAAGAGCTTGGCGGGAATGGGCTGACCGCTTCTCTGTGCTTTACGGTATCGCCGCTCCCGATTGCGAGCGCATCGCCTTC
223▶ A T R D I A E E L G G E W A D R F L V L Y G I A A P D S Q R I A F
7800 TATCGCCTTCTTGACGAGTTCTTCTGA^{EcoRI (7860)}CGGGACTCTGGGGTTCGAAATGACCGACCAAGCGAATTCGCTAGGATTATCCCTAATACCTGCCACCCCACT
257▶ Y R L L D E F F •
7900 CTTAATCAGTGGTGAAGAACGGTCTCAGAAGTCTTTGTTTCAATTGGCCATTTAAGTTTAGTAGTAAAAGACTGGTTAATGATAACAATGCATCGTAAA
^{ScaI (8087)}
8000 ACCTTCAGAAGGAAAGGAGAATGTTTTGTGGACCACTTTGGTTTTCTTTTTTTCGGTGTGGCAGTTTTAAGTTATTAGTTTTTAAAATCAGTACTTTTTAA
8100 TGGAAACAACCTGACCAAAAATTTGTCACAGAATTTTGAGACCCATTAATAAGTTAATGAGAAACCTGTGTGTTCTTTGGTCAACACCGAGACATTT
8200 AGGTGAAAGACATCTAATTCTGTTTTACGAATCTGGAACTTCTGAAAATGTAATCTTGAGTTAACACTTCTGGGTGGAGAATAGGGTTGTTTTCCC
8300 CCCACATAATTGGAAGGGGAAGGAATATCATTAAAGCTATGGGAGGGTTGCTTTGATTACAACACTGGAGAGAAATGCAGCATGTTGCTGATTGCCTGT
8400 CACTAAAACAGGCCAAAACCTGAGTCTTGGGTTGCATAGAAAGCTG
▶