

# pVITRO2-hygro-GFP/SEAP

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

Catalog code: pvitro2-gfpssp

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

For research use only

Version 20H19-MM

## PRODUCT INFORMATION

### Contents

- 20 µg of pVITRO2-hygro-GFP/SEAP provided as lyophilized DNA
- 1 ml Hygromycin B Gold at 100 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 Hygromycin B Gold 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 plasmids developed mainly for *in vitro* studies. They allow the ubiquitous and constitutive co-expression of two genes of interest. pVITRO plasmids can be stably transfected in mammalian cells and the genes of interest are expressed at high levels. Each pVITRO plasmid is available with either two multiple cloning sites or two reporter genes.

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

pVITRO2-hygro-GFP/SEAP also can be used for cloning of open reading frames (ORF). Both reporter genes are flanked by unique sites (BspHI/AvrII for GFP and NcoI/NheI 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.

### Hygromycin B usage:

This antibiotic can be used for *E. coli* at 50-100 µg/ml in liquid or solid media and at 50-500 µg/ml to select Hygromycin-resistant mammalian cells.

## 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 protein<sup>1</sup>. 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<sup>2</sup>. The enhancement varies from 2-fold in non-permissive cells to 20-fold in permissive cells.

- **CMV enhancer:** The major immediate early enhancer of the human cytomegalovirus (HCMV), located between nucleotides -118 and -524, is composed of unique and repeated sequence motifs. The HCMV enhancer can substitute for the 72-bp repeats of SV40 and is severalfold more active than the SV40 enhancer<sup>3</sup>.

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

- **GFP gene:** This red-shifted variant of the jellyfish GFP gene encodes a green fluorescent protein that absorbs blue light (major peak at 480 nm) and emits green light (major peak at 505 nm).

- **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 expression<sup>4</sup>.

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

- **hph gene** confers resistance to Hygromycin B both in *E. coli* and mammalian cells. In bacteria, *hph* is expressed from the constitutive *E. coli* EM7 promoter. In mammalian cells, *hph* is transcribed from the CAG 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.*<sup>5</sup>

1. Eisenstein RS. & Munro HN. 1990. Translational regulation of ferritin synthesis by iron. *Enzyme* 44(1-4):42-58.
2. Dean D.A. *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* 141(2):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(14):2697-700.
5. Carswell S. & Alwine J.C. 1989. Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. *Mol. Cell Biol.* 10:4248-58.

## TECHNICAL SUPPORT

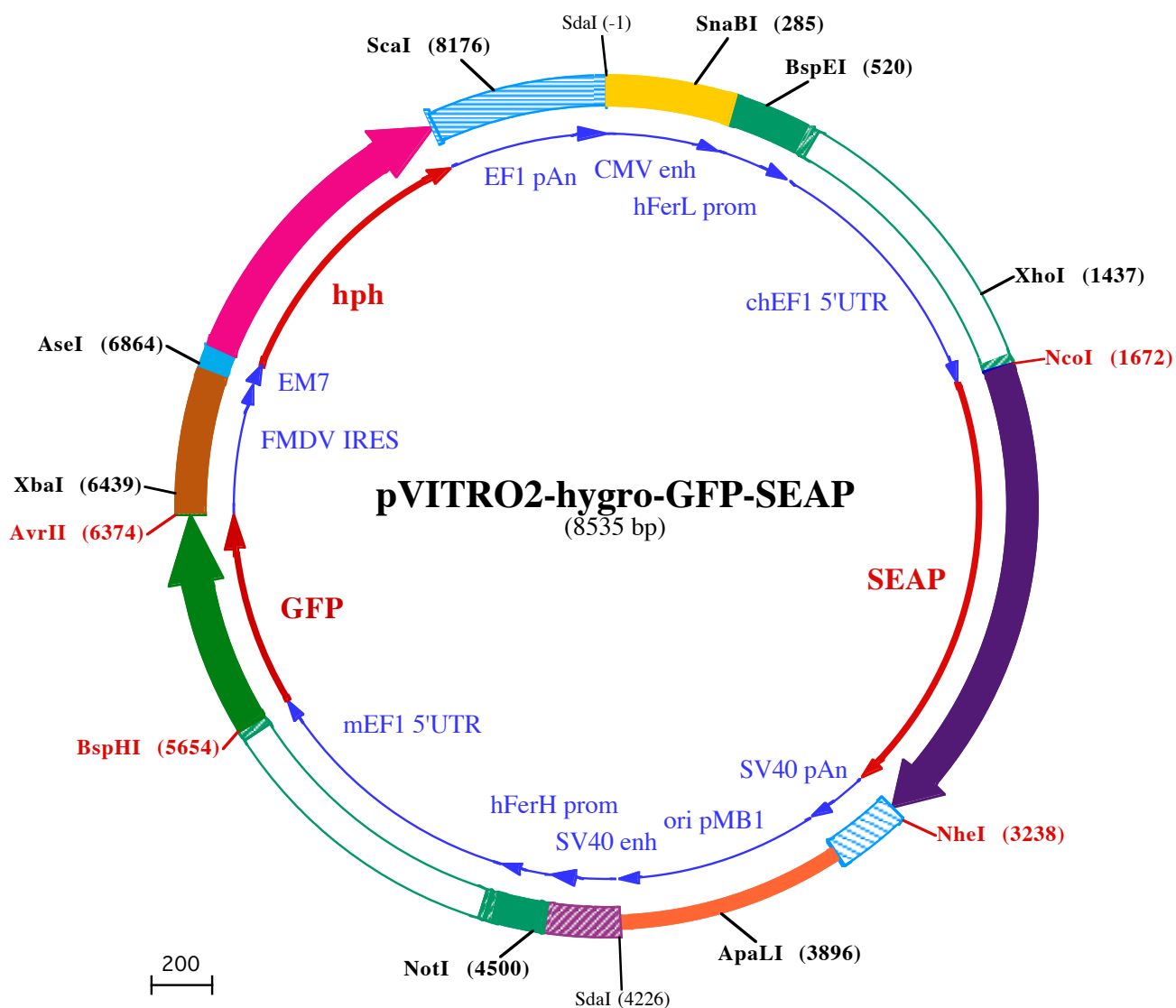
InvivoGen USA (Toll-Free): 888-457-5873

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](mailto:info@invivogen.com)



SdaI (-1)  
1 CCTGCAGGCGTTACATAACTTACGGTAAATGGCCCGCTGGCTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAA

101 CGCCAATAGGGACTTTCCATTGACGTCAATGGTGGAGTATTTACGGTAAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTA

SnaBI (285)  
203 TTGACGTCAATGACGGTAAATGGCCCGCTGGCATTATGCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGCTA

305 TTACCATGATGATGCGGTTTTGGCAGTACATCAATGGCGTGGATAGCGGTTTGACTCACGGGGATTTCGAAGTCTCCACCCCAATTGACGTCAATGGGAGT

406 TTGTTTTGACTAGTCAGGGCCCCAACCCCCCAAGCCCCATTTCACAACACGCTGGCGCTACAGGCGGTGACTTCCCTTGCTTTGGGGGGGGGGCTG

BspEI (520)  
507 AGACTCCTATGTGCTCCGGATTGGTCAGGCACGGCCTTCGGCCCGCTCTGCCACCGCAGATTGGCCGCTAGGCCTCCCGAGCGCCCTGCCTCCGAGG

608 GCGCGGCACCATAAAAGAAGCCGCCCTAGCCACGTCCCTCGCAGTTCGGCGGTCCCGCGGTCTGTCTCAAGCTTGCCGCAGAACACAGGtaagtgc

709 gtgtgtggttcccgcgggcctggcctctttacgggttatggcccttgctgcttgaattacttccatgcccctggctgcagtagctgattcttgatcccca

811 gcttcgggttgaagtgggtgggagagttcgaggccttgcgcttaaggagccccttcgctcgtgcttgagttgagcctggcttggcgctggggcgccg

913 cgtgctaactcgtgtggcaccttcgcgctgtctcgtgctttcgctaagtctctagccatttaaaattttgataaccagctgcgacgcttttttctggcg

1015 agatagctcttgaatatgcccccaagatctgcacactggtatctcggtttttggggcgcgggcgagcgggcccgtgctcccagcgcacatgctcgg

1116 cgaggcggggcctgcgagcggccaccgagaatcggaagggttagtctcaaaactggccggcctgctcgggtgcttggcctcgcgcccgctgtatcgccc

1218 cgccctggggcgcaaggctggcccggcggcaccagttgctgagcggaaagatggccgcttcccggccctgctgcaggagctcaaaatggaggacgcggg

1320 gccgggagagcgggggggtgagtcaccacacaaaaggaaaaggcctttccttctcatccgtcgcttcatgtgactccacggagtagccggcgccgtcca

XhoI (1437)  
1422 ggcacctcgattagttctcgagcttttggagtacgtcgtctttaggttggggggaggggttttatgcatggagtttccccacactgagtggtgggagact

1523 gaagagttaggccagcttggcacttgatgtaattctccttgaatttgcctttttgagtttgatcttgcctcattctcaagcctcagacagtggttcaaa

NcoI (1672)  
1625 gttttttcttccatttccagGTGTCGTGAAAACCTACCCCTAAAAGCCACCATGTTCTGGGGCCTGCATGCTGCTGCTGCTGCTGCTGGCCTGAGGC

1727 TACAGCTCTCCCTGGGCATCATCCAGTTGAGGAGGAGAACCCGACTTCTGGAACCGGAGGCAGCCGAGGCCCTGGGTGCCCAAGAAGCTGCAGCCTG

181▶ 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 A K K L Q P

1829 CACAGACAGCCGCAAGAACCTCATCATCTTCTGGCGATGGGATGGGGGTGTCTACGGTGACAGCTGCCAGGATCCTAAAAGGGCAGAAGAAGGACAAAC

52▶ 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 G Q K K D K

1931 TGGGGCCTGAGATACCCCTGGCTATGGACCGCTCCCATATGTGGCTCTGTCCAAGACATACAATGTAGACAAACATGTGCCAGACAGTGGAGCCACAGCCA

86▶ 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 P D S G A T A

2033 CGGCCTACCTGTGCGGGGTCAAGGGCAACTTCCAGACATTGGCTTGTAGTGCAGCCGCCGCTTAAACCAGTGCAACACGACACCGGCAACGAGGTCTATCT

120▶ 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 T T R G N E V I

2135 CCGTGATGAATCGGGCAAGAAAGCAGGGAAGTCAAGTGGGAGTGGTAACCACCACACGAGTGCAGCAGCCCTCGCCAGCCGGCACCTACGCCACACGGTGA

154▶ 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 A G T Y A H T V

2237 ACCGCAACTGGTACTCGGACGCCGACGTGCCTCGCCCGCCAGGAGGGGTGCCAGGACATCGCTACGAGCTCATCTCCAACATGGACATTGATGTGA

188▶ 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 L I S N M D I D V

2339 TCCTGGTGGAGCCGAAAGTACATGTTTCGCATGGGAACCCGACCCCTGAGTACCCAGATGACTACAGCCAAAGTGGGACCAGGCTGGACGGGAAGAATC

222▶ I L G G G R K Y M F R M G T P D P E Y P D D Y S Q G G T R L D G K N

2441 TGGTGCAGGAATGGCTGGCGAAGCGCCAGGGTCCCGGTATGTGTGGAACCGCACTGAGCTCATGCAGGCTTCCCTGGACCCGCTGTGACCCATCTCATGG

256▶ L V Q E W L A K R Q G A R Y V W N R T E L M Q A S L D P S V T H L M

2543 GTCTCTTTGAGCCTGGAGACATGAAATACGAGATCCACCGAGACTCCACACTGGACCCCTCCCTGATGGAGATGACAGAGGCTGCCCTGGCGCTGCTGAGCA

290▶ G L F E P G D M K Y E I H R D S T L D P S L M E M T E A A L R L L S

2645 GGAACCCCGCGGCTTCTTCTCTTCTGTGGAGGGTGGTGCATCGACCAGGTGCATCACGAAAGCAGGGCTTACCGGGCACTGACTGAGACGATCATGTTCC

324▶ R N P R G F F L F V E G G R I D H G H H E S R A Y R A L T E T I M F

2747 ACGACGCCATTGAGAGGGCGGGCCAGCTACCCAGCGAGGAGGACAGCTGAGCCTGTCACTGCCGACCACTCCACGCTTCTCTTCCGAGGCTACCCCC

358▶ D D A I E R A G Q L T S E E D T L S L V T A D H S H V F S F G G Y P

2849 TGCGAGGAGCTCATCTTCGGCTGGCCCTTGCAAGGCCCGGACAGGAAGCCTACACGGTCTCTATACGAAACGGTCCAGGCTATGTGCTCAAGG

392▶ L R G S S I F G L A P G K A R D R K A Y T V L L Y G N G P G Y V L K

2951 ACGGCGCCCGCGGATGTTACCGAGAGCGAGAGCGGGAGCCCGAGTATCGGACGAGTCAAGTGCAGGCTGACGAGGCTGCCCTGGACGAAGAGACCCACGAGGCGAGGACG

426▶ D G A R P D V T E S E S G S P E Y R Q Q S A V P L D E E T H A G E D

3053 TGGCGGTGTTTCGCGCGGCCCGCAGGCGCACCTGGTTCACGGCGTGCAGGAGCAGACCTTCATAGCGCACGTCATGGCCTTCGCCGCTGCTGGAGCCCT  
460▶ V A V F A R G P Q A H L V H G V Q E Q T F I A H V M A F A A C L E P  
NheI (3238)  
3155 ACACCGCTGCGACCTGGCGCCCCCGCGCACACCACCGACCGCGCACCCGGGGCGGTCCCGGTCCAAGCGTCTGGATTGAAGCTAGCTGGCCAGACATG  
494▶ Y T A C D L A P P A G T T D A A H P G R S R S K R L D •  
3257 ATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATT  
3359 ATAAGCTGCAATAACAAGTTAACAACAACAATTGCATTCAATTTATGTTTCAGGTTTCAGGGGGAGGTGTGGGAGGTTTTTTAAAGCAAGTAAACCTCTAC  
3461 AAATGTGGTATGGAATGTTAATTAAGTACCATGACCAAAATCCCTTAACGTGAGTTTTCTGTTCCACTGAGCGTCAGACCCCGTAGAAAAGATCAAAGGA  
3562 TCTTCTTGAGATCCTTTTTTTCTGCGCGTAATCTGCTGCTTGCAAACAAAAAACACCCTACCAGCGGTGGTTTGTGGCCGATCAAGAGCTACCAACT  
3664 CTTTTCCGAAGGTAAGTGGCTTCAGCAGAGCGCAGATACCAATACTGTTCTTCTAGTGTAGCCGTAGTTAGGCCACCCTTCAAGAAGCTCTGTAGCACCG  
3766 CCTACATACCTCGCTCTGCTAATCCTGTTACCAGTGGCTGCTGCCAGTGGCGATAAGTCGTGTCTTACCGGGTTGGACTCAAGACGATAGTTACCGGATAAG  
ApaLI (3896)  
3868 GCGCAGCGGTGGGGCTGAACGGGGGGTTCGTGCACACAGCCAGCTTGGAGCGAACGACCTACCCGAAGTACCTACAGCGTGAGCTATGAGAAAGC  
3970 GCCACGCTTCCCGAAGGGAGAAAGCGGACAGGTATCCGGTAAGCGGCAGGGTCCGAACAGGAGAGCGCACGAGGGAGCTTCCAGGGGAAACGCCCTGGTAT  
4072 CTTTATAGTCTGTGCGGTTTCGCCACCTCTGACTTGAGCGTCGATTTTTGTGATGCTCGTCAGGGGGCGGAGCCTATGGAAAAACGCCAGAACCGGGCC  
SdaI (4226)  
4174 TTTTTACGGTTCCTGGCCTTTTGTGTCCTTTGCTCACATGTTCTTAATTAACCTGCAGGGCCTGAAATAACCTCTGAAAGAGGAAGTGGTTAGGTA  
4273 CCTTCTGAGGCTGAAAGAACCAGCTGTGGAATGTGTGTCAGTTAGGGTGTGAAAGTCCCAGGCTCCCAGCAGGCAGAGTATGCAAGCATGCATCTC  
4374 AATTAGTCAGCAACCAGGTGTGAAAGTCCCAGGCTCCCAGCAGGCAGAGTATGCAAGCATGCATCTCAATTAGTCAGCAACCATAGTCCCACTAGT  
NotI (4500)  
4475 TCCGCCAGAGCGCGGAGGGCTCCAGCGCCGCCCTCCCCACAGCAGGGGGCGGGTCCCGGCCACCAGGAGCGGGCTCGGGCGGGCGGGCGCTG  
4577 ATTGGCCGGGGCGGGCTGACGCCGACGCGGTATAAGAGACCACAAGCAGCCGACGGCCAGACGTTCTTCGCCGAAGCTTGCCGTCAGAAGCAGGTTG  
4678 AGGGGCGGGTGTGGCTTCGCGGGCGCCGAGCTGGAGTCTGCTCCGAGCGGGCGGGCCCCGCTGTCGTCGGCGGGGATTAGCTGCGAGCATTCCCGCT  
4780 TCGAGTTGCGGGCGGCGGGAGGAGGAGTGCAGGCTAGCGCAACCCGTAAGCCTGCCTCGTGTCCGGCTTAGGGCTAGCGTGGTGTCCGCGCCGCG  
4882 GCCGCGTCTACTCCGGCCGACTCTGGTCTTTTTTTTTTTTGTGTTGTTGCCCTGCTGCCTTCGATTGCCGTTACGCAATAGGGGCTAACAAAGGGAGGG  
4984 TCGGGGCTTGCTCGCCGGAGCCGGAGAGGTATGGTTGGGAGGAATGGAGGGACAGGAGTGGCGGCTGGGGCCCCGCCCTTCGGAGCACATGTCCG  
5086 ACGCCACCTGGATGGGGCAGGCTGGGGTTTTTCCGAAGCAACCAGGCTGGGGTTAGCGTGCCGAGGCCATGTGGCCCCAGCACCCGGCAGCATCTGGCT  
5188 TGGCGCGCCGCTTGCCTGCCTCCCTAACTAGGGTGAAGCCATCCCGTCCGGCACCAGTTGCGTGCGTGAAAGATGGCCGCTCCCGGGCCTGTTGCAA  
5290 GGAGCTCAAAATGGAGGACGCGCAGCCCGTGGAGCGGGCGGGTGAAGTCAACCAAGGAAGAGGGCCTGGTCCCTACCGGCTGCTGCTTCTCTGTGA  
5392 CCCCCTGGTCTATCGCCGCAATAGTCACTCGGGCTTTTGAACAGGCTAGTGCAGCGGGGGAGGGATGTAATGGCGTTGGAGTTTGTTCACATTTG  
5494 GTGGGTGGAGACTAGTCAGGCCAGCCTGGCGCTGGAAGTCAATTTTGAATTTGTCCTTGTAGTTTTGAGCGGAGCTAATTCTCGGGCTTCTTAGCGGTT  
BspHI (5654)  
5596 AAAGGTATCTTTTAAACCCTTTTGTAGTGTGTGAAAACCACCCTAATTCAAAGCAATCATGAGCAAGGGAGAAAGAACTCTTTACTGGTGTGTCCTCAAT  
1▶ M S K G E E L F T G V V P I  
5698 TCTGTTGAGCTGGATGGTGTGTAATGGCCACAAATCTCTGTGTCTGGTGAAGGTGAAGGAGATGCAACTTATGAAAGCTGACTCTGAAGTTCATTTG  
14▶ L V E L D G D V N G H K F S V S G E G E G D A T Y G K L T L K F I C  
5800 TACAACAGGAAAGCTGCCAGTGCCTTGGCCAACCTCTGGTACCACCCTGACTTATGGTGTCAATGTTTTCAGCAGGTACCCTGACCACATGAAGCAGCATGA  
48▶ T T G K L P V P W P T L V T T L T Y G V Q C F S R Y P D H M K Q H D  
5902 CTTCTTTAAATCTGCAATGCCAGAAGTTATGTTTTCAGGAGAGGACAATCTTCTTAAGGATGATGGAATTTATAAGACAAGGGCAGAAGTGAAGTTGAAG  
82▶ F F K S A M P E G Y V Q E R T I F F K D D G N Y K T R A E V K F E G  
6004 TGATACTGGTTAACAGAATTGAGCTGAAAGCATTGATTTTAAAGGAAGTGAACATTCTGGGTACAAGCTGGAGTACAATAATCTCACAAATG  
116▶ D T L V N R I E L K G I D F K E D G N I L G H K L E Y N Y N S H N V  
6106 TTACATTATGGCAGATAAGCAGAAGAATGGAATTAAGTTAATTTCAAGATTAGACACAACATTGAGGATGGATCTGTCCAAGTGGCAGACCATTACCAGCA  
150▶ Y I M A D K Q K N G I K V N F K I R H N I E D G S V Q L A D H Y Q Q  
6208 GAACACCCTATTGGTGTGAGTGGCCAGTCTCCTCCAGATAATCACTATCTCCGCACTCAATCTGCTGTGCAAGACCCCTAATGAGAAAAGAGACCACAT  
184▶ N T P I G D G P V L L P D N H Y L R T Q S A L S K D P N E K R D H M

AvrII (6374)

6310 GGTCTCTCTGGAGTTTGTGACAGCAGCAGGAATTACTCTGGGAATGGATGAGCTGTACAAGTAACTAGGAGCAGGTTTCCCAATGACACAAAACGTGCA

218▶ V L L E F V T A A G I T L G M D E L Y K •

XbaI (6439)

6412 ACTTGAAACTCCGCTGGTCTTTCCAGGTCTAGAGGGTAACACTTTGTACTGCGTTTGGCTCCACGCTCGATCCACTGGCGAGTGTTAGTAACAGCACTGT

6514 TGCTTCGTAGCGGAGCATGACGGCCGTGGAACTCCTCCTTGGTAACAAGACCCACGGGGCCAAAAGCCACGCCACACGGGCCCGTCATGTGTGCAACCC

6616 CAGCACGGCGACTTTACTGCGAAACCCACTTTAAAGTGACATTGAACTGGTACCCACACACTGGTGACAGGCTAAGGATGCCCTTCAGGTACCCCGAGGTA

6718 ACACGCGACACTCGGGATCTGAGAAGGGGACTGGGGCTTCTATAAAAGCGCTCGGTTTAAAAAGCTTCTATGCCTGAATAGGTGACCGGAGGTCGGCACCTT

AseI (6864)

6820 TCCTTTGCAATTACTGACCCTATGAATACAACTGACTGTTTGACAATTAATCATCGGCATAGTATATCGGCATAGTATAATACGACTCACTATAGGAGGG

6920 CCACCATGAAGAACTGAACTGACAGCAACTTCTGTTGAGAAGTTTCTCATTGAAAAATTTGATTCTGTTTCTGATCTCATGCAGCTGTCTGAAGGTGAAG  
1 M K K P E L T A T S V E K F L I E K F D S V S D L M Q L S E G E

7022 AAAGCAGAGCCTTTCTTTTGGATGTTGGAGGAAGAGTTATGTTCTGAGGGTCAATTCCTGTGCTGATGGTTTTTACAAAGACAGATATGTTTACAGACACT

33▶ E S R A F S F D V G G R G Y V L R V N S C A D G F Y K D R Y V Y R H

7124 TTGCCTCTGCTGCTGCTGCAATCCAGAAGTTCTGGACATTGGAGAATTTCTGAATCTCTCACCTACTGCATCAGCAGAAGAGCACAAGGAGTCACTCTCC

67▶ F A S A A L P I P E V L D I G E F S E S L T Y C I S R R A Q G V T L

7226 AGGATCTCCCTGAACTGAGCTGCCAGCTGTTCTGCAACCTGTTGCTGAAGCAATGGATGCCATTGCAGCAGCTGATCTGAGCCAACTCTGGATTTGGTC

101▶ Q D L P E T E L P A V L Q P V A E A M D A I A A A D L S Q T S G F G

7328 CTTTTGGTCCCAAGGCATTGGTCAGTACACCACTTGGAGGGATTTCAATTTGTCCATTGCTGATCCTCATGTCTATCACTGGCAGACTGTGATGGATGACA

135▶ P F G P Q G I G Q Y T T W R D F I C A I A D P H V Y H W Q T V M D D

7430 CAGTTTCTGCTTCTGTTGCTCAGGCACTGGATGAACTCATGCTGTGGGCAGAAGATTGTCTGAAGTCAGACACCTGGTCCATGCTGATTTTGAAGCAACA

169▶ T V S A S V A Q A L D E L M L W A E D C P E V R H L V H A A D F G S N

7532 ATGTTCTGACAGACAATGGCAGAATCACTGCAGTCAATTGACTGGTCTGAAGCCATGTTTGGAGATTCTCAATATGAGGTTGCCAACATTTTTTTTTGGAGAC

203▶ N V L T D N G R I T A V I D W S E A M F G D S Q Y E V A N I F F W R

7634 CTTGGCTGGCTTGCATGGAACAACAACAAGATATTTTAAAAGAAGACCCAGAAGTGGTGGTCCCCAGACTGAGAGCCTACATGCTCAGAATTGGCC

237▶ P W L A C M E Q Q T R Y F E R R H P E L A G S P R L R A Y M L R I G

7736 TGGACCAACTGTATCAATCTCTGGTTGATGAAACTTTGATGATGCTGCTTGGGCACAAGGAAGATGTGATGCCATTGTGAGGTCTGGTCTGAACTGTTG

271▶ L D Q L Y Q S L V D G N F D D A A W A Q G R C D A I V R S G A G T V

7838 GAAGAACTCAAATGCAAGAAGGTCTGCTGCTGTTGGACTGATGGATGTTGAAATCTGGCTGACTCTGAAAACAGGAGACCTCCACAAGACCCAGAG

305▶ G R T Q I A R R S A A V W T D G C V E V L A D S G N R R P S T R P R

7940 CCAAGGAATGAA TATTAGCTAGATTATCCCTAATACCTGCCACCCACTCTTAATCAGTGGTGAAGAACGGTCTCAGAAGTGTGTTTCAATTGGCCA

339▶ A K E •

8040 TTTAAGTTTAGTAGTAAAAGACTGGTTAATGATAACAATGCATCGTAAAACCTTCAGAAGGAAAGGAGAATGTTTTGTGGACCACCTTGGTTTTCTTTTTG

ScaI (8176)

8142 CGTGTGGCAGTTTTAAGTTATTAGTTTTTAAAATCAGTACTTTTTAATGAAACAACCTTGACCAAAAATTTGTCACAGAATTTTGTGACCCATTAAAAAGT

8244 TAAATGAGAAACCTGTGTGTTCCCTTTGGTCAACACCGAGACATTTAGGTGAAAGACATCTAATCTGGTTTTACGAATCTGAAACTTCTTGAAAAATGTAAT

8346 TCTTGAGTTAACTTCTGGGTGAGAATAGGGTTGTTTTCCCCCACATAATTGGAAGGGGAAGGAATATCATTAAAGCTATGGGAGGGTTGCTTTGAT

8447 TACAACACTGGAGAGAAATGCAGCATGTTGCTGATTGCCTGTCACTAAAACAGGCCAAAAACTGAGTCCTTGGTTGCATAGAAAGCTG