

pVITRO2-hygro-GFP/LacZ

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

Catalog code: pvitro2-gfplacz

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

For research use only

Version 19A21-MM

PRODUCT INFORMATION

Contents

- 20 µg of pVITRO2-hygro-GFP/LacZ 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/LacZ contains the GFP and LacZ reporter genes and can be used as a control vector.

pVITRO2-hygro-GFP/LacZ also can be used for cloning of open reading frames (ORF). Both reporter genes are flanked by unique sites (BspH I/Avr II for GFP and Nco I/Nhe I for LacZ) that allow for convenient cloning of ORF's.

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.

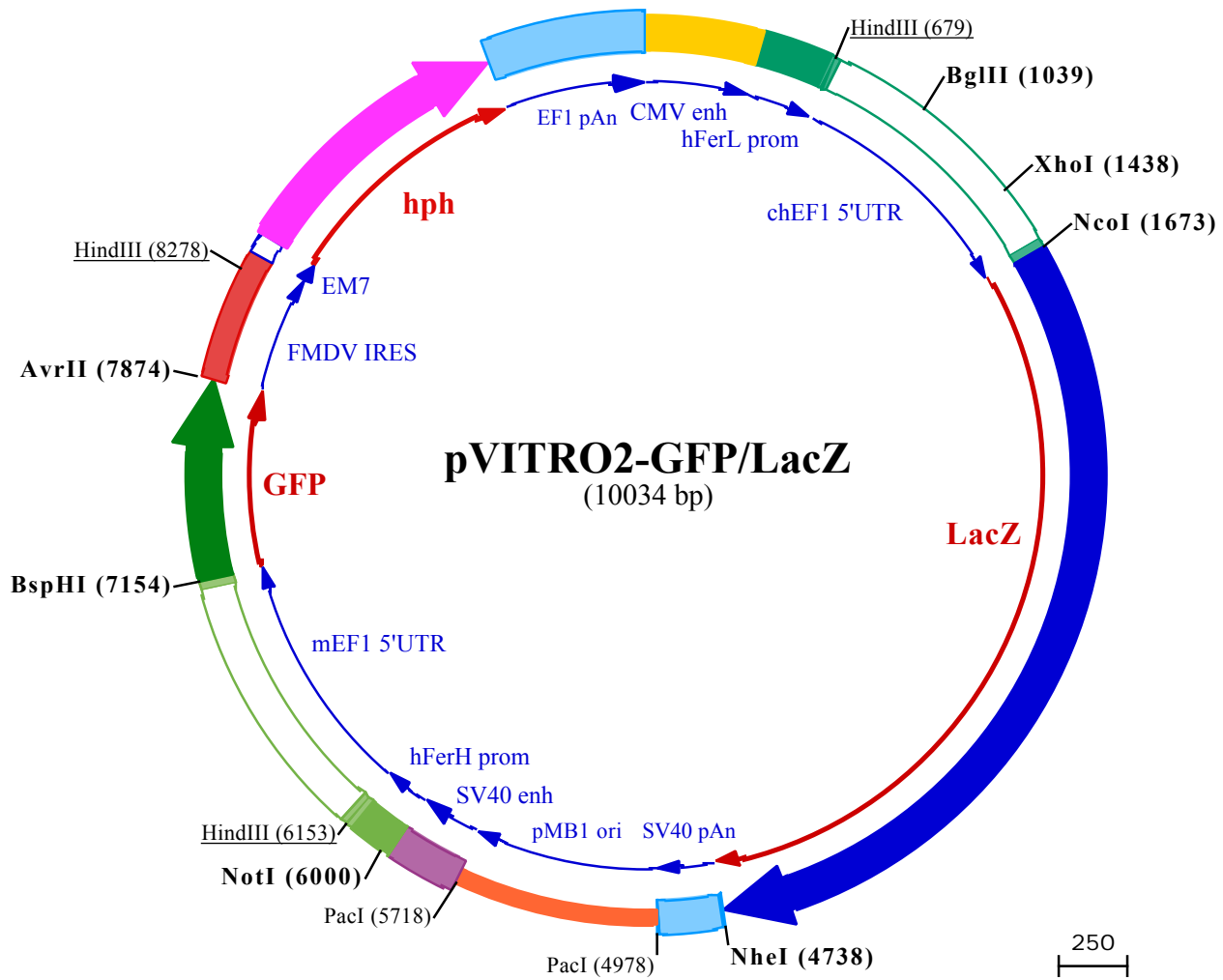
PLASMID FEATURES

- **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 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 range². 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 enhancer³.
- **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 expression⁴.
- **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.
- **LacZ gene:** The *E. coli* lacZ gene codes for the enzyme β-galactosidase which catalyzes the hydrolysis of the substrate X-Gal to produce a blue color that is easily visualized under a microscope.
- **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. 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



1 CCTGCAGGCGTTACATAACTTACGGTAAATGGCCCGCTGGCTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAA
101 CGCCAATAGGGACTTTCCATTGACGTCAATGGGTGGAGTATTTACGGTAAACTGCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCC
201 TATTGACGTCAATGACGGTAAATGGCCCGCTGGCATTATGCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATC
301 GCTATTACCATGATGATGCGGTTTTGGCAGTACATCAATGGGCGTGGATAGCGGTTTGAICTACGGGGATTTCGAAGTCTCCACCCCATTTGACGTCAATG
401 GGAGTTTGTTTGACTAGTCAGGGCCCCAACCCCCCAAGCCCCATTTACAACACGCTGGCGCTACAGGCGGTGACTTCCCCTTGCTTTGGGGCGGG
501 GGGCTGAGACTCCTATGTGCTCCGGATTGGTCAGGCACGGCCTTCGGCCCCGCTCTGCCACCGCAGATTGGCCGCTAGGCCTCCCCGAGCGCCTGCC
601 TCCGAGGGCCGGCCACCATAAAAGAAGCCGCTAGCCAGTCCCCTCGCAGTTCGGCGGTCCCGGGTCTGTCTCAAGCTTCCGCCAGAACACAGg
701 taagtgcctgtgtggttcccggggcctggcctctttacgggttatggccttgcgtgccttgaattacttccatgccctggctgcagtacgtgat t c
801 ttgatcccagcctcgggttgggaagtgggtgggagagttcgaggccttgcgcttaaggagccccttgcctcgtgcttgagttgaggcctggcttgggcg
901 ctggggccgcccgtgctaactctggtggcaccttcgpcctgtctcgctgcttctgctaagtctctagccatttaaaat ttttgataaccagctgcgagc
1001 cttttttctggcgagatagcttctgtaaaggggccaagctgcacactggtat ttcggtttttggggccgggggggcgacggggcccgtgcgtccc
1101 agcgcacatgttcggcgaggcggggcctgcgagcgggcccaccgagaatcggacgggggtagctctcaaatggccgacctgctctggtgcctggcctcgc
1201 gccgacctgtatgccccgcctggggcgcaaggctggcccgttcggcaccagttgcgtgagcggaaagatggccgcttcccggcctgctgcagggagc
1301 tcaaatggaggacgcccggcgcccgggagagcggggcggtgagt caccacacaaaggaaaaggccttctctcctcatccgtcgcttcatgtgactcca
1401 cggagtaccggggcgcgtccaggcacctcgattagttctcgagcttttgagtagctgctcttaggttggggggagggttttatgcatggagtttcc
1501 ccacactgagtggtggagactgaagagttaggcccagcttggcacttgatgtaattctccttggaaatttgccttttgagtttggatcttgctcattc
1601 tcaagcctcagacagtggttcaaagttttttcttccatttcagGTGTCGTGAAACTACCCCTAAAAGCCACCATTGGACCCTGTTGTGCTGCAAAGGAG
1701 AGACTGGGAGAACCCTGGAGTGACCCAGCTCAACAGACTGGCTGCCACCCTCCCTTTGCCTCTTGGAGGAACCTGAGGAAGCCAGGACAGACGGCCC
1801 AGCCAGCAGCTCAGGCTCTCAATGGAGAGTGGAGGTTTGCCTGGTTCCTGCCCTGAAGCTGTGCTGAGTCTGGCTGGAGTGTACCTCCCAGAGG
1901 CTGACACTGTTGTGGTGGCCAGCAACTGGCAGATGCATGGTATGATGCCCCATCTACACCAATGTCACCTACCCATCACTGTGAACCCCTTTTGT
2001 GCCACTGAGAACCCACTGGCTGTACAGCCTGACCTTCAATGTTGATGAGAGTGCTGCAAGAAGGCCAGACCAGGATCATCTTTGATGGAGTCAAC
2101 TCTGCCTTCCACCTCTGGTGCATGGCAGGTGGTGGCTATGGCCAAGACAGCAGGCTGCCCTCTGAGTTGACCTCTCTGCCTTCCAGAGCTGGAG
2201 AGAACAGGCTGGCTGCATGGTGCATGGTCTGATGGCAGCTACCTGGAAGACCAAGCATGTGGAGGATGCTGGCATCTTCAGGGATGTGAGCCT
2301 GCTGCACAAGCCACCACCCAGATTTCTGACTTCCATGTTGCCACCAGGTTCAATGATGACTTCAGCAGAGCTGTGCTGGAGGCTGAGGTGCAGATGCTG
2401 GGAGAATCAGAGACTACCTGAGAGTACAGTGGCCTTGGCAAGTGGAGCCAGGCTGGCCTTGGCAGCCCTTGGAGGAGAGATCATTGATG
2501 AGAGAGGAGGCTATGCTGACAGAGTACCCTGAGGCTCAATGTGGAGAACCCCAAGTGTGGTCTGCTGAGATCCCCAACCTCTACAGGCTGTTGTTGA
2601 GCTGCACACTGCTGATGGCACCCTGATGAAGCTGAAGCCTGTGATGTTGGATTGAGAGAAGTCAGGATTGAGAATGGCCTGCTGCTCAATGGCAAG
2701 CCTCTGCTCATCAGGGGAGTCAACAGGCATGAGCACCACCCTCTGCATGGACAAGTATGGTGAACAGACAATGGTGAAGATATCTGCTAATGAAGC
2801 AGAACAACTTCAATGTCTGCTGCTGCTACTACCCCAACCACCCTCTGCTGACACCCTGTGTGACAGGATGGCCTGTATGTTGTTGATGAAGCCAA
2901 CATTGAGACACATGCCATGGTCCCATGAACAGGCTCACAGATGACCCAGGTTGCCTGCCATGTCTGAGAGAGTACCCAGGATGGTGCAGAGAC
3001 AGGAACACCCCTCTGTGATCATCTGGTCTCTGGCAATGAGTCTGGACATGGAGCAACCATGATGCTCTCTACAGGTGGATCAAGTCTGTTGACCCCA
3101 GCAGACCTGTGCAGTATGAAGGAGTGGAGCAGACCCACAGCAGACATCATCTGCCCATGTATGCCAGGTTGATGAGGACCAGCCCTCCCTGC
3201 TGTGCCAAGTGGAGCATCAAGAAGTGGCTCTCTGCTGGAGAGACCAGCCTCTGATCTGTGTAATGCACATGCAATGGGCAACTCTCTGGGA
3301 GCCTTTGCCAAGTACGGCAAGCTTCAGACAGTACCCAGGCTGCAAGGAGGATTTGTGTGGACTGGTGGACCAATCTCTCATCAAGTATGATGAGA
3401 ATGGCAACCCCTGCTGCTATGGAGGAGACTTTGGTGACACCCCAATGACAGGCAGTTCTGCATGAATGGCTGCTCTTTCAGACAGGACCCTCA
3501 CCCTGCCCTCAGAGGGCCAAAGCACCAGCAACAGTTCTCCAGTTGAGGCTGCTGGACAGACCATTGAGGTGACATCTGAGTACCTTTCAGGCACTCT
609 sProAlaLeuThrGluAlaLysHisGlnGlnGlnPhePheGlnPheArgLeuSerGlyGlnThrIeGluValThrSerGluTyrLeuPheArgHisSer

3601 GACAATGAGCTCCTGCATGGATGGTGGCCCTGGATGGCAAGCCTCTGGCTTCTGGTGAGGTGCTCTGGATGTGGCCCTCAAGGAAAGCAGCTGATTG
643▶ AspAsnGluLeuLeuHisTrpMetValAlaLeuAspGlyLysProLeuAlaSerGlyGluValProLeuAspValAlaProGlnGlyLysGlnLeuIeG
3701 AACTGCCTGAGCTGCCTCAGCCAGAGTCTGTGGACAACTGTGGCTAACAGTGAAGGTTGTCAGCCCAATGCAACAGCTTGGTCTGAGGCAGGCCACAT
676▶ IuLeuProGluLeuProGlnProGluSerAlaGlyGlnLeuTrpLeuThrValArgValValGlnProAsnAlaThrAlaTrpSerGluAlaGlyHisI
3801 CTCTGCATGGCAGCAGTGGAGGCTGGGTGAGAACCCTCTCTGTGACCCTGCCTGCCTCTCATGCCATCCCTCACCTGACAACATCTGAAATGGACTTC
709▶ eSerAlaTrpGlnGlnTrpArgLeuAlaGluAsnLeuSerValThrLeuProAlaAlaSerHisAlaIeProHisLeuThrThrSerGluMetAspPhe
3901 TGCATTGAGCTGGGCAACAAGAGATGGCAGTTCAACAGGCAGTCTGGCTTCTGTCTCAGATGTGGATTGGAGACAAGAAGCAGCTCTCACCCCTCA
743▶ CyslIeGluLeuGlyAsnLysArgTrpGlnPheAsnArgGlnSerGlyPheLeuSerGlnMetTrpIeGlyAspLysLysGlnLeuLeuThrProLeuA
4001 GGGACCAATTACCAGGGCTCCTCTGGACAATGACATTGGAGTGTCTGAGGCCACCAGGATTGACCCAAATGCTTGGGTGGAGAGGTGGAAGGCTGCTGG
776▶ rgAspGlnPheThrArgAlaProLeuAspAsnAspIeGlyValSerGluAlaThrArgIeAspProAsnAlaTrpValGluArgTrpLysAlaAlaGl
4101 ACCTACCAGGCTGAGGCTGCCCTGCTCAGTGCACAGCAGACACCCCTGGCTGATGCTGTTCTGATCACCACAGCCATGCTTGGCAGCACCAAGGCAAG
809▶ yHisTyrGlnAlaGluAlaAlaLeuLeuGlnCysThrAlaAspThrLeuAlaAspAlaValLeuIeThrThrAlaHisAlaTrpGlnHisGlnGlyLys
4201 ACCCTGTTTATCAGCAGAAAGACCTACAGGATTGATGGCTCTGGACAGATGGCAATCACAGTGGATGTGGAGTTGCCCTGACACACCTCACCTGCAA
843▶ ThrLeuPheIeSerArgLysThrTyrArgIeAspGlySerGlyGlnMetAlaIeThrValAspValGluValAlaSerAspThrProHisProAlaA
4301 GGATTGGCCTGAACTGCAACTGGCACAGGTGGCTGAGAGGGTGAAGTGGCTGGGCTTAGGCCCTCAGGAGAATACCCTGACAGGCTGACAGCTGCCTG
876▶ rglIeGlyLeuAsnCysGlnLeuAlaGlnValAlaGluArgValAsnTrpLeuGlyLeuGlyProGlnGluAsnTyrProAspArgLeuThrAlaAlaCy
4401 CTTTGACAGTGGGACCTGCCTCTGTCTGACATGTACCCCTTATGTGTTCCCTTCTGAGAATGGCTGAGGTGTGGCACCAGGGAGCTGAATATGGT
909▶ sPheAspArgTrpAspLeuProLeuSerAspMetTyrThrProTyrValPheProSerGluAsnGlyLeuArgCysGlyThrArgGluLeuAsnTyrGly
4501 CCTCACCAGTGGAGGGGAGACTTCCAGTTCAACATCTCCAGTACTCTCAGAACAGCTCATGGAAACCTCTCACAGGCACCTGCTCCATGCAGAGGAGG
943▶ ProHisGlnTrpArgGlyAspPheGlnPheAsnIeSerArgTyrSerGlnGlnLeuMetGluThrSerHisArgHisLeuLeuHisAlaGluGly
4601 GAACCTGGCTGAACATTGATGGCTTCCACATGGGCATTGGAGGAGATGACTCTTGGTCTCCTCTGTGTCTGCTGAGTTCAGTTATCTGCTGGCAGGTA
976▶ lyThrTrpLeuAsnIeAspGlyPheHisMetGlyIeGlyGlyAspAspSerTrpSerProSerValSerAlaGluPheGlnLeuSerAlaGlyArgTy

NheI (4738)

4701 CCACTATCAGCTGGTGTGGTGCCAGAAGTAAACCTGAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCAACTAGAATGCAG
1009▶ rHisTyrGlnLeuValTrpCysGlnLys•••
4801 TGAAAAAATGCTTTATTGTGAAATTTGTGATGCTATTGCTTTATTGTAAACCATTATAAGCTGCAATAAACAAGTTAAACAACAACAAATTGCATTCAAT

PacI (4978)

4901 TTATGTTTCAGGTTTCAGGGGAGGTGTGGAGGTTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAAATGTTAATTAAGTAGCCATGACCAAA
5001 ATCCCTTAACTGAGTTTTCGTTCCACTGAGCGTCAGACCCCTAGAAAAGATCAAAGGATCTTCTTGAGATCCTTTTTTTCTGCGCGTAATCTGCTGCT
5101 TGCAAAACAAAAAACCCCGTACCAGCGGTGGTTTTGTTGCCGATCAAGAGCTACCAACTCTTTTTCCGAAGGTAAGTGGCTTCAGCAGAGCCGAGAT
5201 ACCAAATACTGTTCTTCTAGTGTAGCCGTAGTTAGGCCACCACTCAAGAACTCTGTAGACCCGCTACATACCTCGCTGCTAATCTGTTACCAGTG
5301 GCTGCTGCCAGTGGCGATAAGTCGTGCTTACCAGGTTGGACTCAAGACGATAGTTACCCGATAAGGCCAGCGGTGGGCTGAACGGGGGGTTTCGTGCA
5401 CACAGCCAGCTTGGAGCGAACGACCTACACCGAACTGAGATACCTACAGCGTGTAGCTATGAGAAAGCGCCAGCTTCCCGAAGGAGAAAGCGGCAG
5501 GTATCCGGTAAGCGGCAGGCTCGAAACAGGAGAGCCAGAGGGAGCTTCCAGGGGAAACGCTGGTATCTTTATAGTCTGTGGGTTTCGCCACCTC
5601 TGACTTGAGCGTCGATTTTGTGATGCTCGTCAGGGGGCGGAGCCTATGAAAAACGCCAGCAACGCGCCTTTTTACGGTTCTTGGCCTTTTGTGCTGC

PacI (5718)

5701 CTTTTGCTCACATGTTCTTAATTAACCTGCAGGCTGAAATAACCTCTGAAAGAGGAACTTGTTAGGTACCTTCTGAGGCTGAAAGAACCAGCTGTGG
5801 AATGTGTGTAGTTAGGTTGGAAAGTCCCAGGCTCCCAGCAGGCAGAAGTATGCAAAGCATGCATCTCAATTAGTCAGCAACCAGGTGTGGAAAGT

NotI (6)

5901 CCCCAGGCTCCCAGCAGGCAGAAGTATGCAAAGCATGCATCTCAATTAGTCAGCAACCATAGTCCCACTAGTTCGCCAGAGCGCGGAGGGCTCCAG
6001 CGGCCGCCCTCCCCACAGCAGGGGCGGGTCCCAGCCACCAGGAGCGGGCTCGGGCGGGCGGCTGATTGGCCGGGGCGGCTGACGCCG

HindIII (6153)

6101 ACGCGCTATAAGAGACCACAAGCGACCCGAGGGCCAGACGTTCTTCGCCGAAGCTTCCCTCAGAACGCAGGTGAGGGCGGGTGTGGCTTCCGCGGG
6201 CCGCCGAGCTGGAGGCTCTGCTCCGAGCGGGCGGGCCCGCTGCTGCGCGGGGATTAGTCTGCGAGCATTCCCGCTTCGAGTTGCGGGCGGGCGGGA
6301 GGCAGAGTGGCAGGCCTAGCGGCAACCCGTAGCCTCGCTCGTGTCCGGCTTGAAGCCTAGCGTGGTGTCCGCGCCGCCGCGCTGCTACTCCGGCCG
6401 CACTCTGGTCTTTTTTTTTTTTGTGTTGTTGTCCTGCTGCCTTCGATTGCCGTTACAGAAAGGGGCTAACAAAGGAGGGTGGCGGGCTTGTCTGCC
6501 GGAGCCCGGAGAGGTCATGTTGGGGAGGAATGGAGGGACAGGAGTGGCGGCTGGGCGCCCGCCCTTCGGAGCACATGTCCGACGCCACCTGGATGGG
6601 GCGAGGCTGGGTTTTTCCGAAAGCAACAGGCTGGGTTAGCGTGCCGAGGCCATGTGGCCCCAGCACCCGGCACGATCTGGCTTGGCGGGCGCGCT
6701 TGCCCTGCCTCCCTAACTAGGGTGGAGCCATCCCGTCCGGCACCAGTTGCGTGGTGGAAAGATGGCCGCTCCCGGGCCCTGTTGCAAGGAGCTCAAAAT
6801 GGAGGACGCGGCAGCCCGGTGGAGCGGGCGGGTGGTACCCACACAAAGAAAGAGGGCCTGGTCCCTCACCAGGCTGCTGCTTCTGTGACCCCGTGGTC
6901 CTATCGGCCGCAATAGTACCTCGGGCTTTTGTAGCAGCGGTAGTCCGGCGGGGGAGGGGATGTAATGGCGTTGGAGTTTGTTCACATTTGGTGGGTGG
7001 AGACTAGTCAGGCCAGCTGGCGCTGGAAGTCATTTTTGGAATTTGTCCCTTGAGTTTGTAGCGGAGCTAATTCTCGGGCTTCTTAGCGGTTCAAGGT

BspHI (7154)

7101 ATCTTTTAAACCTTTTTTAGGTTGTGAAAACCCAGCTAATTCAAAGCAATCATGAGCAAGGGAGAAGAACTTTTACTGGTGTGTTCCCAATTCTG
▶ MetSerLysGlyGluGluLeuPheThrGlyValValProIeLeu

7201 GTTGAGCTGGATGGTGTGATGTGAATGGCCACAAATCTCTGTGTCTGGTGAAGGTGAAGGAGATGCAACTTATGGAAGCTGACTCTGAAGTTCATTGTGA
16▶ ValGluLeuAspGlyAspValAsnGlyHisLysPheSerValSerGlyGluGlyGluGlyAspAlaThrTyrGlyLysLeuThrLeuLysPheI leCysT
7301 CAACAGGAAAAGCTGCCAGTGCCTTGGCCAACCTCTGGTGACCACCCTGACTTATGGTGTTCATATGTTTCAGCAGGTACCCTGACCACATGAAGCAGCATGA
49▶ hrThrGlyLysLeuProValProTrpProThrLeuValThrThrLeuThrTyrGlyValGlnCysPheSerArgTyrProAspHisMetLysGlnHisAs
7401 CTTCTTTAAATCTGCAATGCCAGAAGGTTATGTTTCAGGAGAGGACAATCTTCTTTAAGGATGATGAAAATTATAAGACAAGGGCAGAAGTGAAGTTGAA
82▶ pPhePheLysSerAlaMetProGluGlyTyrValGlnGluArgThrI lePhePheLysAspAspGlyAsnTyrLysThrArgAlaGluValLysPheGlu
7501 GGTGATACACTGGTTAACAGAATTGAGCTGAAAGGCATTGATTTAAGGAAGATGGAACATCTCGGGTCAACAGCTGGAGTACAACATAAATCTCACA
116▶ GlyAspThrLeuValAsnArgI leGluLeuLysGlyI leAspPheLysGluAspGlyAsnI leLeuGlyHisLysLeuGluTyrAsnTyrAsnSerHisA
7601 ATGTTTACATTATGGCAGATAAGCAGAAGAATGGAATTAAGGTTAATTTCAAGATTAGACACAACATTGAGGATGGATCTGTCCAACCTGGCAGACCATTA
149▶ snValTyrI leMetAlaAspLysGlnLysAsnGlyI leLysValAsnPheLysI leArgHisAsnI leGluAspGlySerValGlnLeuAlaAspHisTy
7701 CCAGCAGAACACCCTATTTGGTGTATGGCCAGTTCTCTCCAGATAATCACTATCTCCGCACTCAATCTGCTCTGTCCAAAGCCCTAATGAGAAAAGA
182▶ rGlnGlnAsnThrProl leGlyAspGlyProValLeuLeuProAspAsnHisTyrLeuArgThrGlnSerAlaLeuSerLysAspProAsnGluLysArg
AvrII (7874)
7801 GACCACATGGTCTCTCTGGAGTTTGTGACAGCAGCAGGAATTACTCTGGGAATGGATGAGCTGTACAAGTAAACCTAGGAGCAGGTTTCCCAATGACAC
216▶ AspHisMetValLeuLeuGluPheValThrAlaAlaGlyI leThrLeuGlyMetAspGluLeuTyrLys•••
7901 AAAACGTGCAACTTGAACTCCGCCTGGTCTTCCAGGTCTAGAGGGGTAACACTTTGTACTGCGTTTGGCTCCACGCTCGATCCACTGGCGAGTGTAG
8001 TAACAGCACTGTTGCTTCGTAGCGGAGCATGACGGCGTGGAACTCTCTTGGTAACAAGGACCCACGGGGCCAAAAGCCACGCCACACGGGCCCT
8101 CATGTGTGAACCCAGCACGGCGACTTACTGCGAAACCACTTTAAAGTGACATTGAAACTGGTACCACACACTGGTGACAGGCTAAGGATGCCCTT
HindIII (8278)
8201 CAGGTACCCCGAGGTAACACGGCAGCTCGGGATCTGAGAAGGGGACTGGGGCTTCTATAAAAGCGCTCGGTTTAAAAGCTTCTATGCCTGAATAGGTG
8301 ACCGGAGGTCGGCACCTTTCCTTTGCAATTAAGTCTGACCTATGAATACAACCTGACTGTTTGACAATTAATCATCGGCATAGTATATCGGCATAGTATAATA
8401 CGACTCACTATAGGAGGGCCACCATGAAGAACTGAAGTACAGCAACTTCTGTTGAGAAGTTTCTCATTGAAAAATTTGATTCTGTTTCTGATCTCAT
▶ MetLysLysProGluLeuThrAlaThrSerValGluLysPheLeuI leGluLysPheAspSerValSerAspLeuMe
8501 GCAGCTGTCTGAAGGTGAAGAAAGCAGAGCCTTTCTTTTGGATGTTGGAGGAAGAGGTTATGTTCTGAGGGTCAATCTTGTGCTGATGGTTTTTACAAA
26▶ tGlnLeuSerGluGlyGluGluSerArgAlaPheSerPheAspValGlyGlyArgGlyTyrValLeuArgValAsnSerCysAlaAspGlyPheTyrLys
8601 GACAGATATGTTTACAGACACTTGGCTCTGCTGCTGCCAATTCAGAAAGTTCTGGACATTGGAGAATTTCTGAATCTCTCACCTACTGCATCAGCA
60▶ AspArgTyrValTyrArgHisPheAlaSerAlaAlaLeuProl leProGluValLeuAspI leGlyGluPheSerGluSerLeuThrTyrCysI leSerA
8701 GAAGAGCACAAGGAGTCACTCTCCAGGATCTCCCTGAAACTGAGCTGCCAGCTGTTCTGCAACCTGTTGCTGAAGCAATGGATGCCATTGACGAGCTGA
93▶ rgArgAlaGlnGlyValThrLeuGlnAspLeuProGluThrGluLeuProAlaValLeuGlnProValAlaGluAlaMetAspAlaI leAlaAlaAlaAs
8801 TCTGAGCCAAACCTCTGGATTTGGTCTTTTGGTCCCAAGGCATTGGTCTGACACCACTTGGAGGATTTTCATTTGTGCCATTGCTGATCCTCATGTC
126▶ pLeuSerGlnThrSerGlyPheGlyProPheGlyProGlnGlyI leGlyGlnTyrThrThrTrpArgAspPheI leCysAlaI leAlaAspProHisVal
8901 TATCACTGGCAGACTGTGATGGATGACACAGTTTCTGCTTCTGTTGCTCAGGCCTGGATGAACTCATGCTGTGGGCAGAAGATTGCTCTGAAGTCAGAC
160▶ TyrHisTrpGlnThrValMetAspAspThrValSerAlaSerValAlaGlnAlaLeuAspGluLeuMetLeuTrpAlaGluAspCysProGluValArgH
9001 ACCTGGTCCATGCTGATTTTGGAAAGCAACAATGTTCTGACAGACAATGGCAGAATCACTGCAGTCAATGACTGGTCTGAAGCCATGTTTGGAGATTCTCA
193▶ isLeuValHisAlaAspPheGlySerAsnAsnValLeuThrAspAsnGlyArgI leThrAlaValI leAspTrpSerGluAlaMetPheGlyAspSerGl
9101 ATATGAGGTTGCCAACATTTTGGAGACCTTGGCTGGCTTGCATGGAACAACAACAAGATATTTGAAAGAAGACCCAGAAGCTGGCTGGTTCC
226▶ nTyrGluValAlaAsnI lePhePheTrpArgProTrpLeuAlaCysMetGluGlnGlnThrArgTyrPheGluArgArgHisProGluLeuAlaGlySer
9201 CCCAGACTGAGAGCCTACATGCTCAGAAATGGCCTGGACCAACTGTATCAATCTCTGGTTGATGAAACTTTGATGATGCTGCTTGGGCACAAGGAAGAT
260▶ ProArgLeuArgAlaTyrMetLeuArgI leGlyLeuAspGlnLeuTyrGlnSerLeuValAspGlyAsnPheAspAlaAlaTrpAlaGlnGlyArgC
9301 GTGATGCCATTTGAGGCTGGTCTGGAAGTCTGGAAGAACTCAAATGCAAGAAGGCTGCTGCTGTTGGACTGATGGATGTTTGAAGTTCTGGC
293▶ ysAspAlaI leValArgSerGlyAlaGlyThrValGlyArgThrGlnI leAlaArgArgSerAlaAlaValTrpThrAspGlyCysValGluValLeuAl
9401 TGACTCTGAAACAGGAGACCTCCACAAGCCAGAGCCAAGGAATGAATATTAGCTAGATTATCCCTAATACCTGCCACCCACTCTTAAATCAGTGGT
326▶ aAspSerGlyAsnArgArgProSerThrArgProArgAlaLysGlu•••
9501 GGAAGAACGGTCTCAGAAGCTGTTGTTCAATTGGCCATTTAAGTTAGTAGTAAAAGACTGGTTAATGATAACAATGCATCGTAAAACCTTCAAGAGGA
9601 AAGGAGAATGTTTTGTGGACCACTTTGGTTTTCTTTTTTGGCGTGGCAGTTTTAAGTTATTAGTTTTTAAAATCAGTACTTTTTAATGAAACAACCTG
9701 ACCAAAAATTTGTCACAGAATTTTGGAGCCATTAAGAAAGTTAAATGAGAAACCTGTGTCTCTTTGGTCAACACCGAGACATTTAGGTGAAAGACAT
9801 CTAATCTGGTTTTACGAATCTGAAACTTCTTGAAGTGAATTTCTGAGTTAACACTTCTGGGTGGAGAATAGGGTTGTTTTCCCCACATAATTGG
9901 AAGGGGAAGGAATATCATTAAAGCTATGGGAGGTTGCTTTGATTACAACACTGGAGAGAAATGCAGCATGTTGCTGATTGCCTGTCACTAAAACAGGC
10001 CAAAACTGAGTCTTGGGTTGCATAGAAAGCTG
▶