

# pDRIVE-h $\beta$ -Actin

A plasmid with a native human  $\beta$ -Actin promoter

Catalog # pdrive-hbact

For research use only

Version # 05A17-SV

## PRODUCT INFORMATION

### Content:

- 1 disk of lyophilized GT100 *E. coli* bacteria transformed by a pDRIVE plasmid.
- GT100 genotype is: *F*-, *mcrA*,  $\Delta$ (*mrr-hsdRMS-mcrBC*),  $\emptyset$ 80*lacZ* $\Delta$ M15,  $\Delta$ *lacX74*, *recA1*, *endA1*.
- 4 pouches of *E. coli* Fast-Media® Zeo

### Shipping and storage:

- Products are shipped at room temperature.
- Transformed bacteria should be stored at -20°C. Bacteria are stable up to one year when properly stored.
- Store *E. coli* Fast-Media® Zeo 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.
- Bacteria have been lyophilized, and their viability upon resuspension has been verified.
- Promoter activity has been confirmed by transient transfection of 293 cells as well as other selected cell lines.

## GENERAL PRODUCT USE

pDRIVE is an expression plasmid containing a native or composite promoter of interest. pDRIVE may be used to:

- **Subclone a promoter of interest into another vector.** Unique restriction sites are present at each end of the promoter allowing convenient excision. The 5' sites are *Sda* I, and *Spe* I. *Sda* I is compatible with *Nsi* I and *Pst* I. *Spe* I is compatible with *Avr* II, *Nhe* I and *Xba* I. The 3' restriction site is *Nco* I which includes the ATG start codon, and is compatible with *Bsp*H I and *Bsp*LU11 I.
- **Compare the activity of different promoters** in transient transfection experiments. Each pDRIVE promoter drives the expression of the *LacZ* reporter gene which allows for testing of the promoter's activity in transient transfection experiments. Furthermore, the *LacZ* gene is flanked by unique restriction sites (*Nco* I and *Eco*R I) for easy replacement with a different gene of interest.

## PROMOTER CHARACTERISTICS

Element	Name	Origin	Size bp
Core promoter	$\beta$ -actin	Human	1454
5'UTR	$\beta$ -actin	Human	933
Enhancer	-	-	-

### $\beta$ -Actin promoter

Beta-actin is a highly conserved protein ubiquitously expressed in all eukaryotic cells as a component of the cytoskeleton and as a mediator of internal cell motility. A 1.2-kb fragment of the  $\beta$ -actin 5' flanking region is sufficient for efficient transcription<sup>1</sup>. This 1.2-kb fragment contains a TATA box, a CCAAT box and two highly conserved elements. This fragment was shown to drive high levels of expression of a transgene *in vitro*<sup>2</sup>. InvivoGen provides a larger fragment that includes the 5'UTR of the  $\beta$ -actin gene.

### References:

1. Sugiyama H. *et al.* 1988. Strong transcriptional promoter in the 5' upstream region of the human beta-actin gene. *Gene* 65(1):135-9
2. Muller SR. *et al.* 1990. Efficient transfection and expression of heterologous genes in PC12 cells. *DNA Cell Biol* 9(3):221-9

## PLASMID FEATURES

- **LacZ gene** encodes  $\beta$ -galactosidase an enzyme that catalyzes the hydrolysis of X-Gal, producing a blue precipitate that can be 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.
- **pMB1 Ori** is a minimal *E. coli* origin of replication with the same activity as the longer Ori.
- **EM7** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*.
- **Sh ble** gene confers zeocin resistance therefore allowing the selection of transformed *E. coli* carrying a pDRIVE plasmid.

**Note:** Stable transfection of clones cannot be performed due to the absence of an eukaryotic promoter upstream of the *Sh ble* gene.

## METHODS

### Growth of pDRIVE-transformed bacteria:

Use sterile conditions to do the following:

- 1- Resuspend the lyophilized *E. coli* by adding 1 ml of LB medium in the tube containing the disk. Let sit for 5 minutes. Mix gently by inverting the tube several times.
- 2- Streak bacteria taken from this suspension on a zeocin LB agar plate prepared with the *E. coli* Fast-Media® Zeo agar provided (see below).
- 3- Place the plate in an incubator at 37°C overnight.
- 4- Isolate a single colony and grow the bacteria in TB supplemented with zeocin using the Fast-Media® Zeo liquid provided (see below).
- 5- Extract the pDRIVE plasmid DNA using the method of your choice.

### Selection of bacteria with *E. coli* Fast-Media Zeo:

*E. coli* Fast-Media® Zeo is a **new, fast and convenient** way to prepare liquid and solid media for bacterial culture by using only a microwave. *E. coli* Fast-Media® Zeo is a TB (liquid) or LB (solid) based medium with zeocin, and contains stabilizers.

*E. coli* Fast-Media® Zeo can be ordered separately (catalog code # fas-zn-1, fas-zn-s).

### Method:

- 1- Pour the contents of a 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.

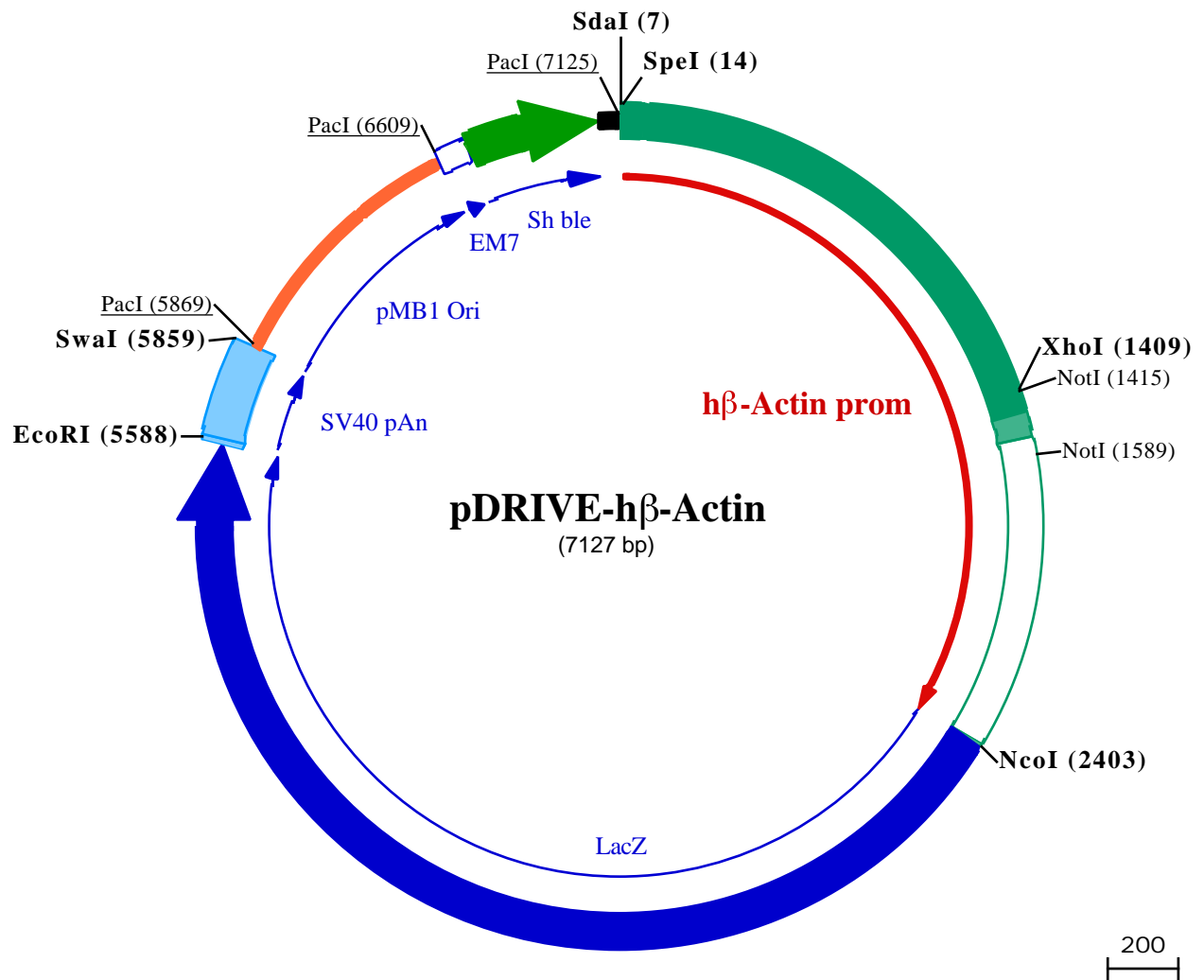
**Note:** Do not reheat solidified Fast-Media® as the antibiotic will be permanently destroyed by the procedure.

## TECHNICAL SUPPORT

Toll free (US): 888-457-5873  
Outside US: (+1) 858-457-5873  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)  
Website: [www.invivogen.com](http://www.invivogen.com)

**InvivoGen**

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



**SdaI (7)**            **SpeI (14)**

1 CCTGCAGGGCCACTAGTTCATGTCTTATATGGACTCATCTTTGCCCTATTGGACACACACTCAATGAACACCTACTACGGCTGCAAAGAGCCCGC  
101 AGGCCTGAGGTGCCCCACCTCACCCTCTTCTATTTTTGTGTA AAAATCCAGCTTCTGTACCACCTCCAAGGAGGGGAGGAGGAGGAAGGCAGGT  
201 TCCTCTAGGCTGAGCGAATGCCCTCTGTGGTCCCACGCCACTGATCGCTGCATGCCACCACCTGGGTACACACAGTCTGTGATTCCCGGAGCAGAAC  
301 GGACCTGCCACCCGGTCTTGTGTCTACTCAGTGGACAGACCCAAGGCAAGAAAGGTGACAAGGACAGGGTCTTCCAGGCTGGCTTTGAGTTCCTA  
401 GCACCGCCCCGCCCAATCTCTGTGGACATGGAGTCTTGGTCCCAGAGTCCCCAGCGGCTCCAGATGGTCTGGGAGGGCAGTTCAGCTGTGGCT  
501 CGCATAGCAGACATAACAAGGACGGTGGGCCAGACCCAGGCTGTGTAGACCCAGCCCCCGCCCGCAGTGCCTAGGTACCCACTAACGCCCCAGG  
601 CCTGGTCTTGGCTGGCGTGACTGTTACCCTCAAAAAGCAGGCAGCTCCAGGGTAAAAGGTGCCCTGCCCTGTAGAGCCACCTTCTTCCAGGGCTGGC  
701 GCTGGGTAGGTTTGTAGCCTTATCACGGGCCACCTCCAGCCACTGGACCCTGGCCCTGCCCTGTCTGGGAGTGTGGTCTCGACTTCTAAGTGG  
801 CCGCAAGCCACTGACTCCCCAACCCACACTCTACCTCTCAAGCCAGGTCTCTCCCTAGTGACCCACCAGCACATTTAGCTAGCTGAGCCCCACAG  
901 CCAGAGTCTCAGGCCCTGCTTTCAGGGCAGTGTCTGAAGTCGCAAGGGGAGTACTGCCTGGCCACTCCATGCCCTCAAGAGCTCCTTCTGCA  
1001 GGAGCGTACAGAACCAGGGCCTGGCACCCGTGCAGACCTGGCCACCCACCTGGGCGCTCAGTGCCAAGAGATGTCCACACCTAGGATGTCCCGC  
1101 GGTGGGTGGGGGCCGAGAGACGGGACGGCCGGGGCAGGCTGGCCATGCGGGGCCAACCAGGACTGCCAGCGTGGGGCCGGGGCCACGGCGC  
1201 GCGCCCCAGCCCCGGGCCAGCACCCCAAGGCGGCAACGCCAAAACCTCTCCCTCTCTCTTCTCAATCTCGCTCTCGCTCTTTTTTTTTTTCGCA  
1301 AAAGGAGGGGAGAGGGGTAAAAAATGCTGCATGTGCGGCAAGCCGGTGTAGTGAGCGGCGGGGCCAATCAGCGTGCCTTCCGAAAGTTCCT

NotI (1415)

**XhoI (1409)**

1401 TTTATGGCTCGAGCGGCCGCGCCCTATAAAAACCCAGCGCGCAGCGCCACCACCGCGAGACCGCTCCGCCCGGAGCACAGAGCTCGC  
1501 CTTTGCCGATCCGCCCGCTCCACACCCGCCGCAgt aagccccggccagccgacggggcaggcggtcacggccccggccgagggcgccggcccc  
1601 ttcgccccgtcgagagcccgcgtctgggcccagcggggggcgcatggggggggaaccggaccgccgtggggggcgcgggagaagcccctgggctccgga  
1701 gatgggggacacccccaccagcttgggagggcgagggccgcgtcgggagggcgctccgggggtgcccgtctcggggcgggggcaaccggcggggtctt  
1801 tgtctgagccgggctcttgccaatggggatcgagggtgggcccggagccccgccaggccccggggctggggcgccattgcgcggtgcgcgctg  
1901 gtcctttgggcgtaactgctgcgcgctgggaattggcgctaattgcgcggtgcgcgctgggactcaaggcgtaactgcgcggtgcttctggggccccg  
2001 ggtgcccgggcctgggctggggcgaaggcgggctcggccggaagggggtggggtcgcccggtcccggggcgcttgccgcaacttctgcccagccgctg  
2101 gccccccgaggggtgtggccgctgctgcgcgcgccgaccggcgctgtttgaaccggggcgagggggctggcggccccggtgggaggggggtggggc  
2201 ctggcttctgcccgcgcccggggagcctccgaccagtgttgccttttatggtaataacggcgccggccggcttcttctgccccaatctgggcg  
2301 cgcgccccgccccctggcgccctaaggactcggcgccggaagtggccagggcgggggcgacctcggtcacagcgcccggtattctcgagCTC

NotI (1589)

**NeoI (2403)**

2401 ACCATGGGGGTTCTCATCATCATCATCATGGTATGGCTAGCATGACTGGTGGACAGAAATGGGTCGGGATCTGTACGACGATGACGATAAGGTAC  
➔ MetGlyGlySerHisHisHisHisHisHisGlyMetAlaSerMetThrGlyGlyGlnGlnMetGlyArgAspLeuTyrAspAspAspLysValP  
2501 CTAAGGATCAGCTGGAGTTGATCCCGTCGTTTACAACGTCGTGACTGGGAAAACCCCTGGCGTTACCCAACCTAATCGCCTTGCAGCACATCCCGCTT  
33➔ roLysAspGlnLeuGlyValAspProValValLeuGlnArgArgAspTrpGluAsnProGlyValThrGlnLeuAsnArgLeuAlaAlaHisProProPh  
2601 CGCCAGCTGGCGTAATAGCGAAGAGGCCCGCACCGATCCGCTTCCCAACAGTTGCGCAGCCTGAATGGCGAATGGCGCTTTGCTGGTTCCGGCACA  
66➔ eAlaSerTrpArgAsnSerGluGluAlaArgThrAspArgProSerGlnGlnLeuArgSerLeuAsnGlyGluTrpArgPheAlaTrpPheProAlaPro  
2701 GAAGCGGTGCCGAAAGCTGGCTGGAGTGGATCTCTCTGAGGCCGATACTGTCTGCTCCCTCAAACCTGGCAGATGCACGGTTACGATGCCCCATCT  
100➔ GluAlaValProGluSerTrpLeuGluCysAspLeuProGluAlaAspThrValValValProSerAsnTrpGlnMetHisGlyTyrAspAlaProIleT  
2801 ACACCAACGTAACCTATCCATTACGGTCAATCCGCGTTTGTTCACGGAGAACCCGAGGGTGTACTCGCTCACATTTAATGTTGATGAAAGCTG  
133➔ yrThrAsnValThrTyrProIleThrValAsnProProPheValProThrGluAsnProThrGlyCysTyrSerLeuThrPheAsnValAspGluSerTr  
2901 GCTACAGGAAGGCAGACGCGAATTATTTTTGATGGCGTTAACTCGCGCTTTCATCTGTGGTGAACGGGCGCTGGGTGGTTACGGCCAGGACAGTCTG  
166➔ pLeuGlnGluGlyGlnThrArgIleIlePheAspGlyValAsnSerAlaPheHisLeuTrpCysAsnGlyArgTrpValGlyTyrGlyGlnAspSerArg  
3001 TTGCCGTCTGAATTTGACCTGAGCGCATTTTACCGCGCCGAGAAAACCGCCTCGCGGTGATGGTGTGCGTTGGAGTGACGGCAGTTATCTGGAAGATC  
200➔ LeuProSerGluPheAspLeuSerAlaPheLeuArgAlaGlyGluAsnArgLeuAlaValMetValLeuArgTrpSerAspGlySerTyrLeuGluAspG  
3101 AGGATATGTGGCGGATGACCGGCATTTCCGTGACGTCTCGTTGCTATAAACCCGACTACACAAATCAGCGATTCCATGTGGCCACTCGCTTAATGA  
233➔ InAspMetTrpArgMetSerGlyIlePheArgAspValSerLeuLeuHisLysProThrThrGlnIleSerAspPheHisValAlaThrArgPheAsnAs  
3201 TGATTTACGCCCGCTGACTGGAGGCTGAAGTTCAGATGTGCGCGAGTTGCGTGACTACCTACGGGTAACAGTTCTTTATGGCAGGTTGAAACGCAG  
266➔ pAspPheSerArgAlaValLeuGluAlaGluValGlnMetCysGlyGluLeuArgAspTyrLeuArgValThrValSerLeuTrpGlnGlyGluThrGln  
3301 AGTCCAGCGCCGCTTTCGGCGGTAATATCGATGAGCGGTGGTTATGCCGATCGCGTACACTACGCTGAACGTCGAAACCCGAAAAC  
300➔ ValAlaSerGlyThrAlaProPheGlyGlyGluIleIleAspGluArgGlyGlyTyrAlaAspArgValThrLeuArgLeuAsnValGluAsnProLysL  
3401 TGTGGAGCGCGAAATCCGAATCTCTATCGTGGGTGGTGAAGTGCACACCGCCGACGGCAGCGTATTGAAGCAGAAGCCTGCGATGTCGGTTCCG  
333➔ euTrpSerAlaGluIleProAsnLeuTyrArgAlaValValGluLeuHisThrAlaAspGlyThrLeuIleGluAlaGluAlaCysAspValGlyPheAr  
3501 CGAGGTGGGATTGAAATGCTGCTGCTGCTGAACGGCAAGCCGTTGCTGATTTCAGGGCTAACCGTACGAGCATCATCTCTGCATGGTCAGGTC  
366➔ gGluValArgIleGluAsnGlyLeuLeuLeuLeuAsnGlyLysProLeuLeuIleArgGlyValAsnArgHisGluHisHisProLeuHisGlyGlnVal

3601 ATGGATGAGCAGACGATGGTGCAGGATATCCTGCTGATGAAGCAGAACAACCTTTAACGCCGTGGCGTGTTCGCATTATCCGAACCATCCGCTGTGGTACA  
400 MetAspGluGlnThrMetValGlnAspIleLeuLeuMetLysGlnAsnAsnPheAsnAlaValArgCysSerHisTyrProAsnHisProLeuTrpTyrT  
3701 CCGTGTGGACCCGCTACGGCCTGTATGTGGTGGATGAAGCAATATTGAAACCCACGGCATGGTGCCAATGAATCGTCTGACCCGATGATCCGGCTGGCT  
433 hrLeuCysAspArgTyrGluLeuTyrValValAspGluAlaAsnIleGluThrHisGlyMetValProMetAsnArgMetValThrAlaThrAspIleIleC  
3801 ACCGGCGATGAGCGAACCGGTAACCGGAATGGTGCAGCGGATCGTAATCACCCGAGTGTGATCATCTGGTCGCTGGGGAATGAATCAGGCCACGGCGCT  
466 uProAlaMetSerGluArgValThrArgMetValGlnArgAspArgAsnHisProSerValIleIleTrpSerLeuGlyAsnGluSerGlyHisGlyAla  
3901 AATCACGACGGCTGTATCGCTGGATCAAATCTGTCTGATCCTTCCCGCCGGTGCAGTATGAAGCGCGGAGCCGACACCAGGCCACCGATATTATTT  
500 AsnHisAspAlaLeuTyrArgTrpIleLysSerValAspProSerHisArgProValGlnTyrGluGlyGlyGlyAlaAspThrThrAlaThrAspIleIleC  
4001 GCCCGATGTACGCGCGCTGGATGAAGACCAGCCCTTCCCGCTGTGCCAAATGGTCCATCAAAAAATGGCTTTCGCTACCTGGAGAGACGCGCCCGCT  
533 ysProMetTyrAlaArgValAspGluAspGlnProPheProAlaValProLysTrpSerIleLysLysTrpLeuSerLeuProGlyGluThrArgProLe  
4101 GATCCTTTGCCAATACGCCACGCGATGGGTAAACAGTCTTGGCGGTTTCGCTAAATACTGGCAGGCGTTCGTCAGTATCCCGCTTACAGGGCGGCTTC  
566 ulLeuLeuGlyGluTyrAlaHisAlaMetGlyAsnSerLeuGlyGlyPheAlaLysTyrTrpGlnAlaPheArgGlnTyrProArgLeuGlnGlyPhe  
4201 GTCTGGGACTGGGTGGATCAGTCTGCTGATTAATATGATGAAAAACGGCAACCCGTGGTGGCTTACGGCGGTGATTTGGCGATACGCCGAACGATCGCC  
600 ValTrpAspTrpValAspGlnSerLeuIleLysTyrAspGluAsnGlyAsnProTrpSerAlaTyrGlyGlyAspPheGlyAspThrProAsnAspArgG  
4301 AGTTCTGTATGAACGGTCTGGCTTTTGGCCAGCCAGCCGCATCCAGCGCTGACGGAAAGCAAAACACCAGCAGCAGTTTTTCCAGTTCGGTTTATCCGG  
633 InPheCysMetAsnGlyLeuValPheAlaAspArgThrProHisProAlaLeuThrGluAlaLysHisGlnGlnGlnPhePheGlnPheArgLeuSerG  
4401 GCAAACCATCGAAGTGACCAGGAAATACCTGTTCGTCATAGCGATAACGAGCTCCTGCAGTGGATGGTGGCGCTGGATGGTAAGCGCTGGCAAGCGGT  
666 yGlnThrIleGluValThrSerGluTyrLeuPheArgHisSerAspAsnGluLeuLeuHisTrpMetValAlaLeuAspGlyLysProLeuAlaSerGly  
4501 GAAGTGCCTCTGGATGTCTGCCACAAGGTAACAGTGTGATGAACTGCCTGAACTACCGCAGCCGGAGAGCGCCGGCAACTCTGGCTCACAGTACGGC  
700 GluValProLeuAspValAlaProGlnGlyLysGlnLeuIleGluLeuProGluLeuProGlnProGluSerAlaGlyGlnLeuTrpLeuThrValArgV  
4601 TAGTGCAACCGAACCGCAGCCATGGTCAGAAGCCGGGCACATCAGCGCTGGCAGCAGTGGCGTCTGGCGAAAACCTCAGTGTGACGCTCCCCCGC  
733 alValGlnProAsnAlaThrAlaTrpSerGluAlaGlyHisIleSerAlaTrpGlnGlnTrpArgLeuAlaGluAsnLeuSerValThrLeuProAlaAl  
4701 GTCCCAGCCATCCCGCATCTGACCACCAGCGAAATGGATTTTTGCATCGAGCTGGTAATAAGCGTTGGCAATTTAACGCCAGTCAAGCTTCTTTTCA  
766 aSerHisAlaIleProHisLeuThrThrSerGluMetAspPheCysIleGluLeuGlyAsnLysArgTrpGlnPheAsnArgGlnSerGlyPheLeuSer  
4801 CAGATGTGGATTGGCGATAAAAAACAACCTGTGACGCCGCTGGCGATCAGTTCACCCGTGCACCGCTGGATAACGACATTTGGCGTAAGTGAAGCGACCC  
800 GlnMetTrpIleGlyAspLysLysGlnLeuLeuThrProLeuArgAspGlnPheThrArgAlaProLeuAspAsnAspIleGlyValSerGluAlaThrA  
4901 GCATTGACCTAACGCCCTGGGTGCAACGCTGGAAGCGCGGGCCATTACCAGGCCAAGCAGCGTGTGTGAGTGCACGGCAGATACACTTGTGTATGC  
833 rglAspProAsnAlaTrpValGluArgTrpLysAlaAlaGlyHisTyrGlnAlaGluAlaAlaLeuLeuGlnCysThrAlaAspThrLeuAlaAspAl  
5001 GGTGCTGATTACGACCGCTCAGCGTGGCAGCATCAGGGGAAAACCTTATTTATCAGCCGAAAACCTACCGGATTGATGTAAGTGGTCAAATGGCGATT  
866 aValLeuIleThrThrAlaHisAlaTrpGlnHisGlnGlyLysThrLeuPheIleSerArgLysThrTyrArgIleAspGlySerGlyGlnMetAlaIle  
5101 ACCGTTGATGTTGAAGTGGCGAGCGATACACCGCATCCGGCGGATTTGGCCTGAACTGCCAGCTGGCGCAGGTAGCAGAGCGGGTAAACTGGCTCGGAT  
900 ThrValAspValGluValAlaSerAspThrProHisProAlaArgIleGlyLeuAsnCysGlnLeuAlaGlnValAlaGluArgValAsnTrpLeuGlyL  
5201 TAGGGCCGCAAGAAAACCTATCCGACCGCCTTACTGCCCTGTTTTGACCGCTGGGATCTGCCATTGTACAGATGTATACCCCGTACGCTTCCCGAG  
933 euGlyProGlnGluAsnTyrProAspArgLeuThrAlaAlaCysPheAspArgTrpAspLeuProLeuSerAspMetTyrThrProTyrValPheProSe  
5301 CGAAAACGGTCTGCGCTGGGACCGCGCAATTGAATATGGCCACACCAAGTGGCGCGGACTTCCAGTTCACACTCAGCCGCTACAGCTCAACAGCAA  
966 rGluAsnGlyLeuArgCysGlyThrArgGluLeuAsnTyrGlyProHisGlnTrpArgGlyAspPheGlnPheAsnIleSerArgTyrSerGlnGlnGln  
5401 CTGATGAAAACCGCATCGCCATCTGCTGCACGGGAAGAAGGCACATGGCTGAATATCGACGTTTCCATATGGGATTTGGTGGCAGCAGCTCTGGA  
1000 LeuMetGluThrSerHisArgHisLeuLeuHisAlaGluGluGlyThrTrpLeuAsnIleAspGlyPheHisMetGlyIleGlyGlyAspAspSerTrpS

EcoRI (5588)

5501 GCCCGTCAGTATCGGCGAATTACAGCTGAGCGCCGGTTCGCTACCATACCAGTTGGTCTGGTGTCAAAAAATAAATCTAGTCCGAGAATTCGCTAGCTC  
1033 erProSerValSerAlaGluLeuGlnLeuSerAlaGlyArgTyrHisTyrGlnLeuValTrpCysGlnLys•••  
5601 GACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTG

5701 TGAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTATTTATGTTTCAGGTTACAGGGG

PacI (5869)

SwaI (5859)

5801 GAGGTGTGGGAGGTTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTAGATCCATTTAAATGTTAATTAAGTACGATGACCAAAATCCCTTAACGTGA

5901 GTTTTCGTTCCACTGAGCGTCAGACCCCGTAGAAAAGATCAAAGGATCTTCTTGAGATCCTTTTTTTCTGCGCGTAATCTGCTGCTTGCAAAACAAAAAA

6001 CCACCGCTACCAGCGGTGGTTTGTGGCGGATCAAGAGCTACCAACTCTTTTTCCGAAGGTAACCTGGCTTACGACAGAGCGCAGATACCAAACTACTGTTCC

6101 TTCTAGTGTAGCCGTAGTTAGGCCACCACTTCAAGAAGCTGTAGCACCAGCTACATACCTCGCTCTGCTAATCCTGTTACCAGTGGCTGCTGCCAGTGG

6201 CGATAAGTCTGTCTTACCGGGTTGGACTCAAGACGATAGTTACCGGATAAGGCCGACGCGTGGGCTGAACGGGGGGTTCGTGCACACAGCCAGCTTG

6301 GAGCGAACACCTACACCGAACTGAGATACCTACAGCGTGAGCTATGAGAAAGCGCCACGCTTCCCGAAGGGAGAAAGCGGACAGGTATCCGGTAAGCC

6401 GCAGGGTCCGAACAGGAGAGCGCAGAGGGAGCTTCCAGGGGAAACGCTGGTATCTTTATAGTCTGTGCGGTTTCGCCACCTCTGACTTGAGCGTGC

6501 ATTTTGTGATGCTCTCAGGGGGCGGAGCCTATGAAAAACGCCAGCAACGCGCCCTTTTTACGGTTCCTGGCCTTTTGTGGCCTTTTGTGCATCATG

PacI (6609)

6601 TCTTAATTAATTTTTCAAAAGTAGTTGACAATTAATCATCGGCATAGTATAATACGACTCACTATAGGAGGGCCATCATGGCCAA

MetAlaLys

6701 GTTGACCACTGCTGCCAGTCTCACAGCCAGGGATGTGGCTGGAGCTGTTGAGTTCTGGACTGACAGGTTGGGGTCTCCAGAGATTTGTGGAGGAT

6801 GACTTTGCAGGTGTGGTTCAGAGATGATGTACCCCTGTTTATCTCAGCAGTCCAGGACCAAGTGGTGCCTGACAACACCCCTGGCTTGGGTGTGGGTGAGAG

6901 GACTGGATGAGCTGTATGCTGAGTGGAGTGGTCTCCACCACTTCCAGGATGCCAGTGGCCCTGCCATGACAGAGATTTGGAGAGCAGCCCTGGGG

7001 GAGAGACTTTGCCCTGAGAGACCCAGCAGCAACTGTGTGCACTTTGTGGCAGAGGAGCAGGACTGAGGATAAGAATGAGTTTCAGAAAAGGGGGCCTG

103 yArgGluPheAlaLeuArgAspProAlaGlyAsnCysValHisPheValAlaGluGlnAsp•••

PacI (7125)

7101 AGTGGCCCTTTTTCAACTTAATTA