

pSELECT-blasti-LacZ

A LacZ-expression plasmid selectable with blasticidin

Catalog code: psetb-lacz

<https://www.invivogen.com/pselect-blasti>

For research use only

Version 20117-MM

PRODUCT INFORMATION

Contents

- 20 µg of pSELECT-blasti-LacZ plasmid 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

pSELECT plasmids are specifically designed for strong and constitutive expression of a gene of interest in a wide variety of cell lines. They allow the selection of stable transfectants and offer a variety of selectable markers. pSELECT plasmids contain two expression cassettes: the first drives the expression of the gene of interest and the second drives the expression of a large choice of dominant selectable markers for both *E. coli* and mammalian cells. They are both terminating with a strong polyadenylation signal (polyA) that separates the two expression cassettes thus preventing any transcription interference. The late SV40 polyA terminates the transcription of the gene of interest while the human β -globin polyA terminates the transcription of the selectable marker.

pSELECT-LacZ plasmids can be used as control vectors or for cloning of an open reading frame, as the LacZ gene is flanked by two unique restriction sites: Nco I at the 5' end that encompasses the Start codon, and Nhe I at the 3' end.

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 DH5 α .

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

First expression cassette

- **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1 α (EF-1 α) core promoter¹ and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat². The EF-1 α promoter exhibits a strong activity and yields long lasting expression of a transgene *in vivo*. The R-U5' has been coupled to the EF-1 α core promoter to enhance stability of RNA.

- **LacZ**: 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 mRNA³.

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

Second expression cassette

- **CMV enh/prom**: The human cytomegalovirus immediate-early gene 1 promoter/enhancer was originally isolated from the Towne strain and was found to be stronger than any other viral promoters.

- **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*. The *bsr* gene is driven by the CMV enhancer/promoter in tandem with the bacterial EM7 promoter allowing selection in both mammalian cells and *E. coli*.

- **β Glo pAn**: The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription⁴.

References

1. Kim D.W. *et al.*, 1990. Use of the human elongation factor 1 alpha promoter as a versatile and efficient expression system. *Gene* 2: 217-223.
2. Takebe Y. *et al.*, 1988. SR alpha promoter: an efficient and versatile mammalian cDNA expression system composed of the simian virus 40 early promoter and the R-U5 segment of human T-cell leukemia virus type 1 long terminal repeat. *Mol. Cell Biol.* 1: 466-472.
3. 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: 4248-4258.
4. Yu J. & Russell J.E., 2001. Structural and functional analysis of an mRNP complex that mediates the high stability of human beta-globin mRNA. *Mol Cell Biol.* 21(17):5879-88.

TECHNICAL SUPPORT

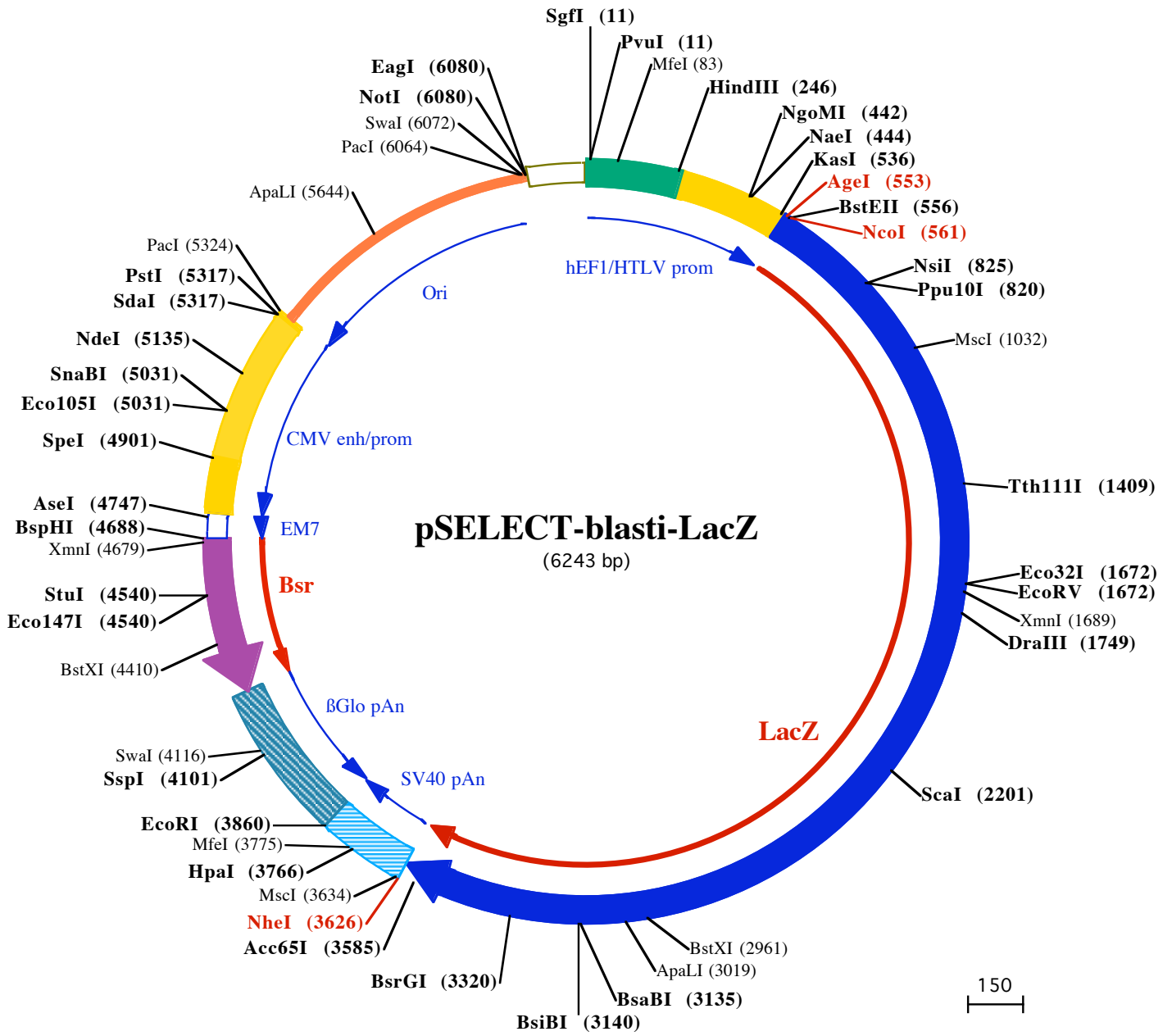
InvivoGen USA (Toll-Free): 888-457-5873

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



PvuI (11) SgfI (11) MfeI (83)

1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGCGAGAGCCACATCGCCACAGTCCCGAGAAAGTTGGGGGAGGGTTCGCAATTGAACGGGTGCCTA

101 GAGAAGTGGCGGGGTAACCTGGAAAGTGATGCTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (246)

201 GTGAACGTTCTTTTCGCAACGGGTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACGCGCCCGCCCTACTGAGGCC

301 GCCATCCACGCGGTTGAGTCGGGTTCTGCCGCCCTCCCGCTGTGGTGCCTCTGAACTGCGTCCGCGCTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMI (442) NaeI (444)

401 GGGCCTTTGTCGGCGCTCCCTTGGAGCTACCTAGACTCAGCGGGCTCTCCACGCTTTGCTGACCTGCTTGTCTCAACTCTACGCTTTTGTCTGTT

NcoI (561) BstEII (556) KasI (536) AgeI (553)

501 TCTGTTCTGGCGGTTACAGATCCAAGCTGTGACCGGGCTACTCTGAGATCACCGGTACCCTGGACCCTGTGTGCTGCAAAAGGAGAGCTGGGAGAA

601 CCCTGGAGTGACCCAGCTCAACAGACTGGCTGCCACCTCCCTTTGCTCTGGAGGAAGCTGAGGAAGCCAGGACAGACAGCCAGCCAGCAGCTC

130 nProGlyVal Thr Gl nLeuAsnArgLeuAl aAl aHi sP roP roPheAl aSer TrpArgAsnSer Gl uGl uAl aArgThrAspArgP roSer Gl nGl nLeu

701 AGGTCCTCAATGGAGAGTGGAGGTTGCTGGTCCCTGCCCTGAAAGCTGTGCCTGAGTCTTGGCTGGAGTGTGACCTCCAGAGGCTGACACTGTTG

470 ArgSer LeuAsnGly uTrpArgPheAl aTrpPheProAl aProGly uAl aVal P roGly uSer TrpLeuGly uCysAspLeuP roGly uAl aAspThr Val V

Ppu10I (820) NsiI (825)

801 TGGTCCAGCACTGGCAGATGCATGGCTATGATGCCCATCTACACCAATGCACCTACCCCATCACTGTGAACCCCTTTTGTGCCACTGAGAA

800 aI Val P roSerAsnTrpGly nMetHisGly TyrAspAl aP roI eTyrThrAsnVal Thr TyrP roI eThr Val AsnP roP roPheVal P roThr Gl uAs

901 CCCCACTGGCTGCTACAGCTGACCTTCAATGTTGATGAGAGCTGGCTGCAAGAGGCCAGACAGGATCATCTTTGATGGAGTCAACTCTGCTTCCAC

1130 nP roThr Gl yCysTyrSer LeuThr PheAsnVal AspGly uSer TrpLeuGly nGl uGly uGl nThr ArgI l eI l ePheAspGly yVal AsnSer Al aPheHis

MscI (1032)

1001 CTCTGGTCAATGGCAGGTGGTGGCTATGGCCAAGCAGCAGGCTGCCCTGAGTTGACCTCTCTGCTTCCCTCAGAGCTGGAGAGAAGCAGGCTGG

1470 LeuTrpCysAsnGly yArgTrpVal Gl yTyrGly uGl nAspSerArgLeuP roSer Gl uPheAspLeuSer Al aPheLeuArgAl aGly yGl uAsnArgLeuA

1101 CTGTCATGGTCTCAGTGGTCTGATGGCAGCTACCTGGAAGACCAAGACATGTGGAGGATGTCTGGCATCTTCAGGGATGTGAGCCTGCTGCACAAGCC

1800 l aVal MetVal LeuArgTrpSerAspGly ySer TyrLeuGly uAspGly nAspMetTrpArgMetSer Gl yI l ePheArgAspVal Ser LeuLeuHisLysP r

1201 CACCACCAGATTTCTGACTTCCATGTTGCCACCAGGTTCAATGATGACTTCAGCAGAGCTGTGCTGGAGGCTGAGGTGCAGATGTGGAGAAGCTCAGA

2130 oThr Thr Gl nI l eSerAspPheHisVal Al aThr ArgPheAsnAspAspPheSer ArgAl aVal LeuGly uAl aGly uVal Gl nMetCysGly yGl uLeuArg

1301 GACTACCTGAGAGTCACAGTGGCCTCTGGCAAGGTGAGACCCAGTGGCTCTGGCACAGCCCTTTGGAGGAGAGATCATTGATGAGAGAGGAGGCT

2470 AspTyrLeuArgVal Thr Val Ser LeuT rpGly nGly uThr Gl nVal Al aSer Gl yThr Al aP roPheGly yGly uI l eI l eAspGly uArgGly yGly T

Tth111I (1409)

1401 ATGCTGACAGAGTCAACCTGAGGCTCAATGTGGAGAAGCCCAAGCTGTGGTCTGCTGAGATCCCAACCTCTACAGGGCTGTGTGGAGCTGCACACTGC

2800 yAl aAspArgVal Thr LeuArgLeuAsnVal Gl uAsnP roLysLeuT rpSer Al aGly uI l eP roAsnLeuTyrArgAl aVal Val Gl uLeuHis sThr Al

1501 TGTGGCAAGCTGATGAAAGCTGAAGCCTGTGATGTGGATTGGATTGAGAAAGTCAAGGATGAGAAATGGCTGCTGCTGCTGCTGCTGCTGCTGCTGCTC

3130 aAspGly yThr LeuI l eGly uAl aGly uAl aCysAspVal Gl yPheArgGly uVal ArgI l eGly uAsnGly yLeuLeuLeuLeuAsnGly yLysP roLeuLeuI l e

EcoRV (1672) Eco32I (1672) XmnI (1689)

1601 AGGGGAGTCAACAGGCATGAGCACCACCTCTGCATGGACAAGTGGATGAACAGACAATGGTGAAGATATCCCTGTAATGAAGCAGAAACAACCTCA

3470 ArgGly yVal AsnArgHis sGly uHi sHi sP roLeuHis sGly yGl nVal MetAspGly uGl nThr MetVal Gl nAspI l eLeuLeuMetLysGly nAsnAsnPheA

DraIII (1749)

1701 ATGCTGTCAAGTGTCTCACTACCCCAACCCCTCTCTGGTACACCTGTGTGACAGGTATGGCCTGTATGTTGTGTGATGAAGCCAACATTGAGACACA

3800 snAl aVal ArgCysSer His sTyrP roAsnHis sP roLeuT rpTyrThr LeuCysAspArgTyrGly yLeuTyrVal Val AspGly uAl aAsnI l eGly uThrHi

1801 TGGCATGGTCCCATGAACAGGCTCACAGATGACCCAGGCTGGCTGCCATGCTGAGAGAGTGAAGGATGACAGGATGGTGAAGCAAGCAACCC

4130 sGly yMetVal P roMetAsnArgLeuThr AspAspP roArgTrpLeuP roAl aMetSer Gl uArgVal Thr ArgMetVal Gl nArgAspArgAsnHis sP ro

1901 TCTGTGATCATCTGGTCTCTGGCAATGAGTCTGGACATGGAGCCAACATGATGCTCTCTACAGTGGATCAAGTCTGTTGACCCAGCAGACCTGTGC

4470 Ser Val I l eI l eT rpSer LeuGly yAsnGly uSer Gl yHi sGly yAl aAsnHis sAspAl aLeuTyrArgTrpI l eLysSer Val AspP roSer ArgP roVal G

2001 AGTATGAAGGAGTGGAGCAGACACCACGCCACAGACATCATCTGCCCATGTATGCCAGGTTGATGAGGACCAGCCCTTCCCTGCTGTGCCAAGTG

4800 l nTyrGly uGly yGly yAl aAspThr Thr Al aThrAspI l eI l eCysP roMetTyrAl aArgVal AspGly uAspGly nP roPheP roAl aVal P roLysTr

ScaI (2201)

2101 GAGCATCAAGAAGTGGCTCTCTGCTGGAGAGACCAGACCTCTGATCCTGTGTAATGCACATGCAATGGGCAACTCTCTGGAGGCTTTGCCAAG

5130 pSer I l eLysLysTrpLeuSer LeuP roGly uThr ArgP roLeuI l eLeuCysGly uTyrAl aHi sAl aMetGly yAsnSer LeuGly yGly yPheAl aLys

2201 TACTGGCAAGCCTTCAAGCAGTACCCAGGCTCAAAGGAGGATTTGCTGGAGCTGGGTGGACCAATCTCTCATCAAGTATGATGAGAATGCAACCCCT

5470 TyrTrpGly nAl aPheArgGly nTyrP roArgLeuGly nGly yGly yPheVal TrpAspTrpVal AspGly nSer LeuI l eLysTyrAspGly uAsnGly yAsnP roT

2301 GGTCTGCATGGAGGAGACTTTGGTGGACACCCCAATGACAGGAGTCTGATGATGATGGCTGGTCTTTGACAGACAGGACCCCTCACCTGCCCTCAC

5800 rpSer Al aTyrGly yGly yAspPheGly yAspThr P roAsnAspArgGly nPheCysMetAsnGly yLeuVal PheAl aAspArgThr P roHi sP roAl aLeuTh

2401 AGAGGCCAAGCACCAGCAAGTCTTCCAGTTCAGGCTGTCTGGACAGACATTGAGGTGACATCTGAGTACCTCTTACGGCACTCTGACAAAGAGCTC

6130 r Gl uAl aLysHis sGly nGly nPhePheGly nPheArgLeuSer Gl yGly nThr I l eGly uVal Thr Ser Gl uTyrLeuPheArgHis sSerAspAsnGly uLeu

2501 CTGCACTGGATGGTGGCCCTGGATGGCAAGCCTCTGGCTTCTGGTGGAGTGGCTCTGGATGGCCCTCAAGGAAGCAGCTGATTGAACCTGCTGAGC

6470 LeuHis sTrpMetVal Al aLeuAspGly yLysP roLeuAl aSer Gl yGly uVal P roLeuAspVal Al aP roGly nGly yLysGly nLeuI l eGly uLeuP roGly uL

2601 TGCCTCAGCCAGAGTCTGCTGGACAAGTGGCTAACAGTGAAGGTTGCTCAGCCCAATGCAACAGCTTGGTCTGAGGAGCCACATCTCTGCATGGCA

6800 euP roGly nP roGly uSer Al aGly yGly nLeuTrpLeuThr Val ArgVal Val Gl nP roAsnAl aThr Al aTrpSer Gl uAl aGly yHi sI l eSer Al aTrpGly

2701 GCACTGGAGGCTGGTGAAGCCTCTCTGACCCCTGGCTGCTGCTCATGCCATCCCTCAGTCAACATCTGAAATGGACTTCTGATGAGCTG

7130 nGly nTrpArgLeuAl aGly uAsnLeuSer Val Thr LeuP roAl aAl aSer His sAl aI l eP roHi sLeuThr Thr Ser Gl uMetAspPheCysI l eGly uLeu

2801 GGCAACAGAGATGGCAGTTCACAGGCAGCTGGCTTCTGCTCAGATGGATTGGAGACAAGAAGCAGCTCTCACCCCTCTCAGGACCAATTC

7470 Gl yAsnLysArgT rpGly nPheAsnArgGly nSer Gl yPheLeuSer Gl nMetTrpI l eGly yAspLysLysGly nLeuLeuThr P roLeuArgAspGly nPheT

BstXI (2961)

2901 CCAGGCTCCTTGGACAATGACATTTGAGTGTCTGAGGCCACAGGATGACCAATGCTTGGTGGAGAGGTGGAAGCTGCTGGACACTACAGGC

7800 hr ArgAl aP roLeuAspAsnAspI l eGly yVal Ser Gl uAl aThr ArgI l eAspP roAsnAl aTrpVal Gl uArgT rpLysAl aAl aGly yHi sTyrGly nAl

ApaLI (3019)

3001 TGAGGCTGCCCTGCTCAGTGCACAGACACCCTGGCTGATGCTGTTCTGATCACCACAGCCATGCTTGGCAGCACAAGCAAGACCTGTTTCATC

8130 aGly uAl aAl aLeuLeuGly nCysThr Al aAspThr LeuAl aAspAl aVal LeuI l eThr Thr Al aHi sAl aTrpGly nHi sGly nGly yLysThr LeuPheI l e

BsiBI (3140) BsaBI (3135)

3101 AGCAGAAAGACCTACAGATTGATGGCTCGACAGATGGCAATCAGATGGATGGAGGTTGCCTCTGACACACCTCACCTGCAAGGATTGGCCTGA

8470 SerArgLysThr TyrArgI l eAspGly ySer Gl yGly nMetAl aI l eThr Val AspVal Gl uVal Al aSer AspThr P roHi sP roAl aArgI l eGly yLeuA

3201 ACTGTCAACTGGCAGAGGTGGCTGAGAGGGTGAACCTGGCTGGGCTTAGGCCCTCAGGAGAATACCCTGACAGGCTGACAGCTGCCTGCTTTGACAGGTG
880▶ snCysGlnLeuAlaGlnValAlaGluArgValAsnTrpLeuGluLeuGlyProGlnGluAsnTyrProAspArgLeuThrAlaAlaCysPheAspArgTrp
BsrGI (3320)

3301 GGACCTGCCTCTGCTGACATGTACACCCCTTATGTGTTCCCTTCTGAGAATGGCCTGAGGTGTGGCACCAGGGAGCTGAACATGGTCTCACCAGTGG
913▶ pAspLeuProLeuSerAspMetTyrThrProTyrValPheProSerGluAsnGlyLeuArgCysGlyThrArgGluLeuAsnTyrGlyProHisGlnTrp
3401 AGGGGAGACTTCCAGTTCAACATCTCCAGGTACTCTCAGCAACAGCTCATGGAAACCTCTCACAGGCACCTGCTCCATGCAGAGAGGGAACTGGCTGA
947▶ ArgGlyAspPheGlnPheAsnIleSerArgTyrSerGlnGlnGlnLeuMetGluThrSerHisArgHisLeuLeuHisAlaGluGluGlyThrTrpLeuA
Acc65I (3585)

3501 ACATTGATGGCTTCCACATGGGCATTGGAGGAGATGACTCTTGGTCTCCTTCTGTGTCTGCTGAGTTCCAGTTATCTGTGGCAGGTACCACATCAGCT
980▶ snIleAspGlyPheHisMetGlyIleGlyGlyAspAspSerTrpSerProSerValSerAlaGluPheGlnLeuSerAlaGlyArgTyrHisTyrGlnLe
MseI (3634)

3601 GGTGTGGTCCAGAAGTAAACCTGAGCTAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGC
1013▶ uValTrpCysGlnLys●●●

NheI (3626)

3701 TTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGTGCATAAACAAGTTAACACAACAATTCATTCTTTTATGTTTCAGG
HpaI (3766) MfeI (3775)

3801 TTCAGGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATCTAAAATACAGCATAGCAAACTTTAACCTCAAAT
EcoRI (3860)

3901 CAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTCA
SspI (4101)

4001 TGGAGTTTAAGATATAGTGTATTTTCCCAAGTTTGAAGTACTAGCTCTTCATTCTTTATGTTTTAAATGCACTGACCTCCCACATTCCCTTTTATGATAAAA

Swal (4116)

4101 TATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTA

4201 GTAGTTGGACTTAGGGAACAAGAAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTAGCTTTAGTTCCTGGTGTACTTGAGGGGGATGAGTT
141▶●●●AsnArgThrTyrLysLeuProIleLeuGlu
4301 CCTCAATGGTGGTTTTGACCAGCTTGCATTATCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTGCATGCCACAGGGGCT
130▶uGluIleThrThrLysValLeuLysGlyAsnMetGluIleLeuValPheCysAspProAlaTyrAspSerIleLeuGluArgCysMetGlyCysProSer
BstXI (4410)

4401 GACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCCTTCTGCCGTTGCTCACAGCAGACCAATG
97▶ValValArgIleSerArgAspValGluAspSerTyrProHisArgValAlaValIleThrAspPheAspLysGlnGlyAsnSerValAlaSerGlyIleA
StuI (4540)

4501 GCAATGGCTTCCAGCACAGACAGTACCCTGCCAATGTAGGCCCTCAATGTGGACAGCAGAGATGATCTCCCAGTCTTGGTCTGTGGCCGCCCGACAT
63▶IleAlaGluAlaCysValThrValArgGlyIleTyrAlaGluIleHisValAlaSerIleIleGluGlyThrLysThrArgIleAlaAlaGlyValHis
Eco147I (4540)

4601 GGTGCTTGTTCCTCATAGAGCATGGTATCTTCTCAGTGGCGACCTCCACAGCTCCAGATCTGCTGAGAGATGTTGAAGGCTTTCATGATGGCCCT
30▶sHisLysAsnAspGluTyrLeuMetThrIleLysGluThrAlaValGluValLeuGluLeuAspGlnGlnSerIleAsnPheThrLysMet
XmnI (4679)

4701 CCTATAGTGAGTCGTATTATACTATGCCGATATACATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCACTAA
AseI (4747)

4801 ACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTGTGATT
SpeI (4901)

4901 ACTAGTCAAAACAACCTCCATTGACGTCAATGGGTGGAGACTTGGAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTAAGTCCAAAACC

SnaBI (5031)

Eco105I (5031)

5001 GCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTT

NdeI (5135)

5101 ACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGCAGTTTACCCTAAATACTCCACCCATTGACGTCAA
5201 TGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTACGCCAGGCGGGCCATTACCCTAA

PaeI (5324)

PstI (5317)

SdaI (5317)

5301 GTTATGTAAACGCCCTGCAGGTTAATTAAGAACATGTGAGCAAAAGCCAGCAAAAGCCAGGAACCGTAAAAAGCCGCGTTGCTGGCGTTTTTCCATAGG
5401 CTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGCGTTTTCCCTCGGAAGCT
5501 CCCTCGTGCCTCTCTGTTCCGACCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTCGGGAAGCGTGCCGCTTCTCATAGCTCACGCTGTAG

ApaLI (5644)

5601 GTATCTCAGTTCGGTGTAGGTGCTTCCGCTCAAGCTGGGCTGTGTGCAGAAACCCCGTTCAGCCGACCGCTGCGCCTTATCCGGTAATATCGTCTT
5701 GAGTCAAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGA
5801 AGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATC
5901 CGGCAAAACAACACCGCTGGTAGCGGTGTTTTTTTTGTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTT

EagI (6080)

6001 ACGGGTCTGACGCTCAGTGAACGAAAACCTCAGCTTAAGGATTTTGGTATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTT
PaeI (6064) Swal (6072) NotI (6080)

6101 ATTTTTCATTACATCTGTGTGTTGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAACGAAACAACAACTAGCAAAATAGGCT
6201 GTCCCCAGTCAAGTGCAGTGCCAGAACATTTCTCTATCGAA