

pBROAD2-LacZ

An optimized vector for mouse and rat transgenesis

Catalog # pbroad2-lacz

For research use only

Version 02B28-MT

PRODUCT INFORMATION

Content:

- 20 µg of pBROAD2-LacZ provided as lyophilized DNA

Storage and Stability:

- Products are shipped at room temperature.
- Lyophilized DNA should be resuspended upon receipt and stored at -20°C (see Methods). Lyophilized DNA is stable 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

The pBROAD2-lacZ plasmid is designed for the expression of *E. coli* β-galactosidase gene in virtually all tissues of transgenic mice and rats. This feature is brought by the ROSA promoter. The murine ROSA26 promoter was initially identified by random retroviral gene trapping in mouse embryonic stem cells¹. This high CpG content promoter was shown to drive ubiquitous expression of the human placental alkaline phosphatase and enhanced green fluorescent protein during embryonic and postnatal development in mouse and rat². The ROSA promoter cloned into pBROAD2-LacZ is the human counterpart of the murine Rosa 26 promoter.

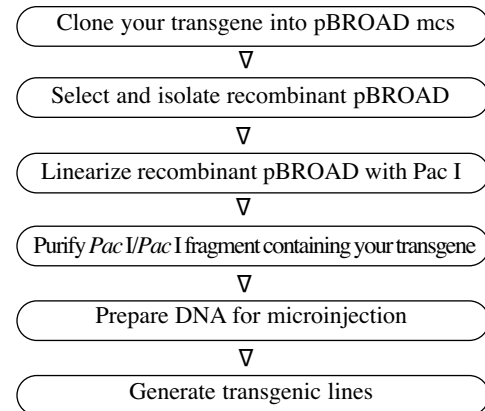
pBROAD2-LacZ expresses a new chemically synthesized gene (*LacZΔCpGNLS*) deprived of CpGs to eliminate interferences of CpG methylation on gene expression. Furthermore, the *E. coli* region is flanked on either side by the well cutting 8 bp-recognizing restriction enzyme *Pac I* that enables linearization and easy excision of the *E. coli* region.

PLASMID FEATURES

- **hROSA prom:** This TATA-less promoter, highly homologous to the murine promoter, was found to be very effective *in vitro* in a very broad range of mammalian cell lines. The strength of the human ROSA promoter is ascribed to the 10 potential Sp1 sites found within the CpG island extending from the core promoter to the first half of 5' untranslated region (5'UTR), the higher number of Sp1 sites never recorded in any natural promoter. The 5'UTR contains an engineered intron of 1200 bp which increases the transcription of the transgene³.
- **LacZ-ΔCpG NLS (pBROAD2-LacZnls):** 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. A nuclear localization signal of SV40 large T has been inserted in the 5' end of the *lacZ* gene to allow the targeting of the chimeric protein to the nucleus. To reduce the immunogenicity of this bacterial gene, InvivoGen has engineered a synthetic *lacZnls* gene that is entirely free of CpG motifs, whereas the wild type *lacZ* gene contains 298 CpG dinucleotides.

- **βGlob pAn:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription⁴.
- **pMB1 ori:** a minimal *E. coli* origin of replication to limit vector size but with the same activity as the longer Ori.
- **Amp:** The ampicillin resistance gene allows the selection of transformed *E. coli* carrying a pBROAD plasmid.

EXPERIMENTAL OUTLINE



METHODS

Plasmid resuspension:

Quickly spin the tube containing the lyophilized plasmid to pellet the DNA. To obtain a plasmid solution at 1µg/µl, resuspend the DNA in 20µl of sterile H₂O. Store resuspended plasmid at -20°C.

Pac I linearization of recombinant pBROAD:

1- Digest 10 µg recombinant pBROAD2 plasmid with 1 to 5 units of *Pac I* restriction enzyme.

Note: *Pac I* may be purchased from New England Biolabs and used at 0.1-0.5 unit per µg plasmid DNA.

2- Incubate at 37°C for 1-2 hours.

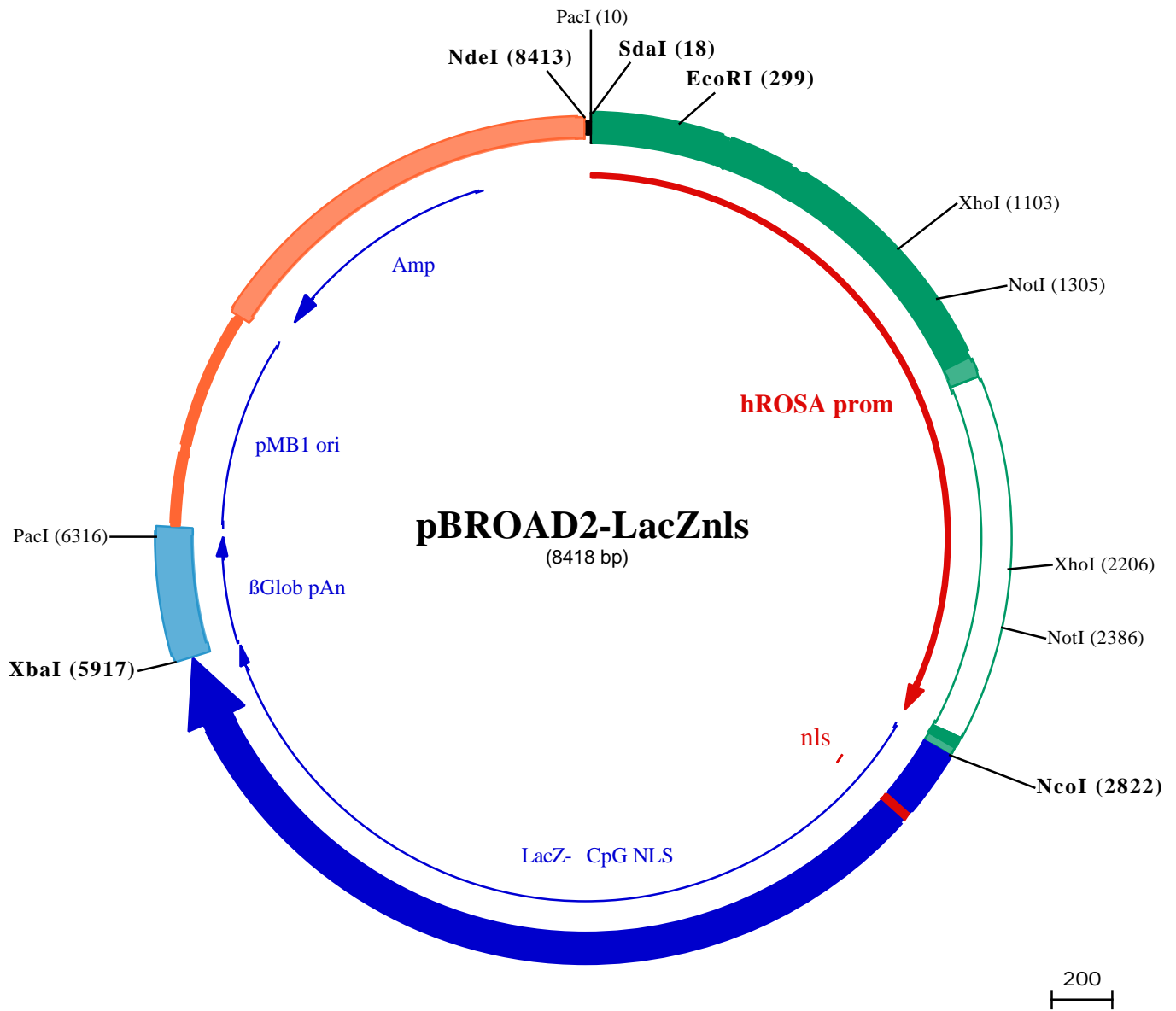
3- Purify the fragment containing the ROSA26 prom-transgene-βGlo pAn cassette by agarose gel following your usual protocol.

References:

1. Zambrowicz BP, et al. 1997. Disruption of overlapping transcripts in the ROSA beta geo 26 gene trap strain leads to widespread expression of beta-galactosidase in mouse embryos and hematopoietic cells. *Proc Natl Acad Sci USA*. 94:3789-94.
2. Kisseberth WC, et al. 1999. Ubiquitous expression of marker transgenes in mice and rats. *Dev Biol*. 214:128-38.
3. Brinster RL, et al. 1988. Introns increase transcriptional efficiency in transgenic mice. *Proc Natl Acad Sci USA* 85(3):836-40
4. Yu J, Russell JE. 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 3-622-34-80
E-mail: info@invivogen.com



PacI (10) SdaI (18)

1 GATCTCGACTTAATTAACCTGCAGGTGAATCATTGCACAAGTAACATGAGAAAGCAGAAAATGCAGGTCATACACGCACCCCTGACCCAGACCAGCAGAG
101 CTGACTGCAGCATCCATATCCAAGAGAAAGACCCTGACGCCCAAGAAGTGAGACAAGCAAGGACTCTATAGAATCAATTAGCATAGAAGGGCTTTCC

EcoRI (299)

201 AACAGTTTAACTTTCCCTCTCATGCGATTTACCTACTTGAACCAGGGCTCTTTCCTACACTCCTCTTACATTCCCGACTTACACGCAGAGGGAAAGAGA
301 ATTCATAAAGGGAATATTTTTCTGCCTTTGAAGATATTCTCACAAGATCGTTCTCCACGCCAAGGCAAGTAAACGACACAATCTGGCTCACTCCAGG
401 CTCGAACCTTACACATTCAACGAGGCTATCTCAGACACGCTGTGGCACACGCCACGGGGAGCCAGAAAACGTGTGGTGGGGGTGGCGAAGGTAATGCCTT
501 TGGGAAGCAGCCATCTGAGGTGGGAGCCAGAAAACGAGAGGGAAGCGTCCAGGAAGATTACGGAGGGGAGATCGCGGCCCCAGAGCGATCAGAGTTG
601 TCTGTCACAAGGCCGCGAGAACCGGGGTAGGGAGTGGGGATCGGGGAGAGAAAAAAGTATGCCTGTGTATTTTCGAGCGGAGGCAGCAAGAGGCCTGT
701 CCTCATTGGAAAGTAAACGTGGAGTAGGCAGTCCCAGGAAAGGGGTGAAGAGGCCTTGGGGAGGGGAAGCGTCTGACCCAGGAAAAACATGAAAG
801 GGGGAGTTGGTTCGCTAGATTAGAGGGGATCTCTCTCCCTGGGAAAATGGGGTGTGCAACGGTGTGTGCAAGCGGCGAGGGGGTGAAGTGGCA
901 GCATCCTCTAAGAGCTTGGGGAGGGCCAGGCCACGCCAAGGAGAGCGAGCGGGGAGACGGAGGAGGTGACCCCTCCCTCCCTCGGGCCCGATC
1001 GTGAGGTTCCGTCTTTTTCTGTCGGACCTTACCTTGTCCAGGCGTGCCTGGGCTGGGCCCGGGTGCAGCGCACGGCACTCCCGGAGGCGCGAG

XhoI (1103)

1101 GACTCGAGTTAGGCCAACCGCGCCACGGCGTTTCTGGCCGGGAATGGCCGTACCCGTGAGGTGGGGGTGGGGGCGAGAAAGCGGAGCGAGCCAA
1201 AGGCGGGGAGGGGGGCGAGGGCCAGGGAAGAGGGGGCCGGCCTACTGTGTTGGCGGACTGGCGGGACTGGGGCTGCGTGTGAGTCTCTGAGCGCAGGGC

NotI (1305)

1301 GGCGGCGCGCCCTCCCCGCGAGCGGGCGGGCGGGCGGGCGGGCGGGCGGGCAGCTCACTAGCCCGCTGCCGAGCGGAAACGCCACTGACCGC
1401 ACGGGGATCCAGCGCGGCGCCAGGGCACCCGGACACGCCCTCCCGCCGCGCATTGGCTCTCCGCCACCGCCCGCACCCATTGGCCACT
1501 CGCCGCAATCAGCGGAAGCCGCGGGCCGCTAGAGAAGAGGCTGTGCTTCCGCGCTCCGGCTCCTCAGAGAGCCTCGGCTAGgt aggggagcggaa
1601 tctggtagggagggagggtgcggtgcactgggggatgggtagggggccgtctggtaggcttgcgggggttccttcccgtgggaagt cgggaaca
1701 taatgtttgttacgttgggagggaaaggggtggctggatgcaggcgggaggaggccccctgcggcaaccggagggggaggagaagggagcggaaaa
1801 tgctcgaaaccggacggagccattgctctcgagagggaggagcgttccggctagcctcttgcgagcttggccgttctcctcccggcgtgtgtgaa
1901 aacacaaatggcgtattctggttggagtaaaactcctgtcagttacaccgtcgggagtagcgagccgttagcgactctcggttgccccctgggtgggg
2001 cgggttaggtaggtaggggtgtagagatgctgggtgtgcggcgcggcccgctcctgcggcgggagggagggtcagtgaattggctctggcgcgggct
2101 cctcccaccctccccttcttccggggagtcggttaccgcggcctgcttgccttcgacacctgattggctgtcgaagctgtgggaccggcccttgc

XhoI (2206)

2201 actggctcaggtctcacatgagcgaaccactgcgcgggagcgggggtggcgggaggcggcgttggtacggtcctccccgagggcggagcgcgcgagt

NotI (2386)

2301 gtctggccccgcccctgcgcaactggcaggaagcgcgctggaggcgggggagggtgcccggccgagacttctggatggcggcgccggcgtccg
2401 cccccgggttcccaccgctgaaggcgagacaagcccagcctgctacaggcactcgtgggggtgggggaggagcgggggtcggtccggctggttgtggg
2501 tgggagcgcctgttctccaaaaaccgagcgagctgcaatcctgaggagctgcggtggaggaggtggagagaagccaccccttctgggaggggga
2601 ggggagtagcgcgaataccttatgggagttcttctgctgcctcccgtcttgaaggaccgctgggctggaagaagccctccctccttctcctcgcg
2701 tgatctcgtcatcgccctccatgctcagctcgttctcgattatgggaggatcttttgcctagacaattgtactaaccttcttctcttctcctcctgac

NcoI (2822)

2801 agGTGTGAAACAGGAAGAGAACCATGGACCCTGTGTGCTGCAAAGGAGAGACTGGGAGAACCTGGAGTGACCAGCTCAACAGACTGGCTGCCACCC
1▶MetAspProValValLeuGlnArgArgAspTrpGluAsnProGlyValThrGlnLeuAsnArgLeuAlaAlaHisPr
2901 TCCCTTTGCCTCTTGGAGGAAGCTGAGGAAGCCAGGACAGAGGCCACCCAGCAGCTCAGTCTCTCAATGGAGAGTGGAGTTTGCCTGGTCCCT
26▶oProPheAlaSerTrpArgAsnSerGluGluAlaArgThrAspArgProSerGlnGlnLeuArgSerLeuAsnGlyGluTrpArgPheAlaTrpPhePro
3001 GCCCTGAAGCTGTGCTGAGTCTTGGCTGGAGTGTGACCTCCAGAGCAGTCCCAAGAAAGAGGAAAGTTGAGGCTGACACTGTGTGGTGCCAA
60▶AlaProGluAlaValProGluSerTrpLeuGluCysAspLeuProGluAlaValProLysLysLysArgLysValGluAlaAspThrValValValProS
3101 GCACTGGCAGATGCATGGCTATGATGCCCCATCTACACCAATGTCACCTACCCATCACTGTGAACCCCTTTTGTGCCACTGAGAACCCCACTGG
93▶erAsnTrpGlnMetHisGlyTyrAspAlaProlIeTyrThrAsnValThrTyrProIeThrValAsnProProPheValProThrGluAsnProThrGl
3201 CTGCTACAGCTGACCTTCAATGTGATGAGAGCTGGCTGCAAGAAGGCCAGACCAGGATCATCTTTGATGGAGTCACTCTGCCTTCCACCTCTGGTG
126▶yCysTyrSerLeuThrPheAsnValAspGluSerTrpLeuGlnGluGlyGlnThrArgIleIePheAspGlyValAsnSerAlaPheHisLeuTrpCys

3301 AATGGCAGGTGGTGGCTATGGCCAAGACAGCAGGCTGCCCTCTGAGTTTGACCTCTCTGCCTTCTCAGAGCTGGAGAGAACAGGCTGGCTGTCATGG
160▶ AsnGlyArgTrpValGlyTyrGlyGlnAspSerArgLeuProSerGluPheAspLeuSerAlaPheLeuArgAlaGlyGluAsnArgLeuAlaValMetV
3401 TGCTCAGGTGGTCTGATGGCAGCTACCTGGAAGACCAAGACATGTGGAGGATGTCTGGCATCTTCAGGGATGTGAGCCTGCTGCACAAGCCCACCACCA
193▶ alLeuArgTrpSerAspGlySerTyrLeuGluAspGlnAspMetTrpArgMetSerGlyl lePheArgAspValSerLeuLeuHisLysProThrThrGl
3501 GATTTCGACTTCCATGTTGCCACCAGGTTCAATGATGACTTCAGCAGAGCTGTGCTGGAGGCTGAGGTGCAGATGTGTGGAGAACTCAGAGACTACCTG
226▶ nl leSerAspPheHisValAlaThrArgPheAsnAspPheSerArgAlaValLeuAlaGluValGlnMetCysGlyGluLeuArgAspTyrLeu
3601 AGAGTCACAGTGGCCCTCTGGCAAGGTGAGACCCAGGTGGCCTCTGGCACAGCCCTTTGGAGGAGAGATCATTGATGAGAGAGGAGGCTATGCTGCACA
260▶ ArgValThrValSerLeuTrpGlnGlyGluThrGlnValAlaSerGlyThrAlaProPheGlyGlyGluI leI leAspGluArgGlyGlyTyrAlaAspA
3701 GAGTCACCCTGAGGCTCAATGTGGAGAACCCCAAGCTGTGGTCTGCTGAGATCCCAACCTCTACAGGGCTGTTGTGGAGCTGCACACTGCTGATGGCAC
293▶ rgValThrLeuArgLeuAsnValGluAsnProLysLeuTrpSerAlaGluI leProAsnLeuTyrArgAlaValValGluLeuHisThrAlaAspGlyTh
3801 CCTGATTGAAGCTGAAGCCTGTGATGTTGGATTTCAGAGAAGTCAGGATTGAGAATGGCCTGCTGCTGCTCAATGGCAAGCCTCTGCTCATCAGGGGAGCT
326▶ rLeul leGluAlaGluAlaCysAspValGlyPheArgGluValArgI leGluAsnGlyLeuLeuLeuAsnGlyLysProLeuLeuI leArgGlyVal
3901 AACAGGCATGAGCACCACCCTCTGCATGGACAAGTGTGGTGAACAGACAATGGTGAAGATATCCTGCTAATGAAGCAGAACAACCTCAATGCTGTCA
360▶ AsnArgHisGluHisHisProLeuHisGlyGlnValMetAspGluGlnThrMetValGlnAspI leLeuLeuMetLysGlnAsnAsnPheAsnAlaValA
4001 GGTGCTCTACTACCCCAACCACCCTCTCTGGTACACCCTGTGTGACAGGTATGGCCTGTATGTTGTTGATGAAGCCAACATTGAGACACATGGCATGGT
393▶ rgCysSerHisTyrProAsnHisProLeuTyrTrpTyrThrLeuCysAspArgTyrGlyLeuTyrValValAspGluAlaAsnI leGluThrHisGlyMetVa
4101 GCCCATGAACAGGCTACAGATGACCCAGGTGGCTGCCATGTCTGAGAGAGTGACCAGGATGGTGCAGAGAGACAGGAACCCCTCTGTGATC
426▶ lProMetAsnArgLeuThrAspAspProArgTrpLeuProAlaMetSerGluArgValThrArgMetValGlnArgAspArgAsnHisProSerVal I le
4201 ATCTGGTCTCTGGCAATGAGTCTGGACATGGAGCAACCATGATGCTCTCTACAGGTGGATCAAGTCTGTTGACCCAGCAGAGCTGTGCAGTATGAAG
460▶ l leTrpSerLeuGlyAsnGluSerGlyHisGlyAlaAsnHisAspAlaLeuTyrArgTrpI leLysSerValAspProSerArgProValGlnTyrGluG
4301 GAGGTGGAGCAGACACCACAGCCACAGACATCATCTGCCCATGTATGCCAGGGTGTATGAGGACCAGCCCTCCCTGCTGTGCCAAGTGGAGCATCAA
493▶ lyGlyGlyAlaAspThrAlaThrAspI leI leCysProMetTyrAlaArgValAspGluAspGlnProPheProAlaValProLysTrpSerI leLy
4401 GAAGTGGCTCTCTGCTGGAGAGACCAGACCTCTGATCCTGTGTGAATATGCACATGCAATGGGCAACTCTCTGGGAGGCTTTGCCAAGTACTGGCAA
526▶ sLysTrpLeuSerLeuProGlyGluThrArgProLeuI leLeuCysGluTyrAlaHisAlaMetGlyAsnSerLeuGlyGlyPheAlaLysTyrTrpGln
4501 GCCTCAGACAGTACCCAGGCTGCAAGGAGGATTTGTGTGGGACTGGGTGGACCAATCTCTCATCAAGTATGATGAGAATGGCAACCCCTGGTCTGCCT
560▶ AlaPheArgGlnTyrProArgLeuGlnGlyPheValTrpAspTrpValAspGlnSerLeuI leLysTyrAspGluAsnGlyLeuProTrpSerAlaT
4601 ATGGAGGAGCTTTGGTGACCCCAATGACAGGCACTCTGCATGAATGGCCTGGTCTTTGACAGCAGGACCCTCAGCCCTCACAGGGCCAA
593▶ yrGlyGlyAspPheGlyAspThrProAsnAspArgGlnPheCysMetAsnGlyLeuValPheAlaAspArgThrProHisProAlaLeuThrGluAlaLy
4701 GCACCAGCAACAGTCTTCCAGTTCAGGCTGTCTGGACAGACCATTGAGGTGACATCTGAGTACCTCTTCAGGCACTCTGACAATGAGCTCCTGCACTGG
626▶ sHisGlnGlnGlnPhePheGlnPheArgLeuSerGlyGlnThrI leGluValThrSerGluTyrLeuPheArgHisSerAspAsnGluLeuLeuHisTrp
4801 ATGGTGGCCTGGATGGCAAGCCTCTGGCTTCTGGTGGAGTGGCCTCTGGATGTGGCCCTCAAGGAAAGCAGCTGATTGAACCTGCCTGAGCTGCCTCAGC
660▶ MetValAlaLeuAspGlyLysProLeuAlaSerGlyGluValProLeuAspValAlaProLeuAspGlnProPheProAlaValProLysTrpSerI leLy
4901 CAGAGTCTGCTGGACAACCTGTGGCTAACAGTGGGTTAGCCCAATGCAACAGCTTGGTCTGAGGAGCCACATCTCTGCATGGCAGCAGTGGAG
693▶ roGluSerAlaGlyGlnLeuTrpLeuThrValArgValValGlnProAsnAlaThrAlaTrpSerGluAlaGlyHisI leSerAlaTrpGlnGlnTrpAr
5001 GCTGGCTGAGAACCCTCTCTGTGACCTGCCTGCTGCCTCTCATGCCATCCCTCACCTGACAACATCTGAAATGGACTTCTGCATTGAGCTGGGCAACAAG
726▶ gLeuAlaGluAsnLeuValThrLeuProAlaAlaSerHisAlaI leProHisLeuThrThrSerGluMetAspPheCysI leGluLeuGlyAsnLys
5101 AGATGGCAGTTCAACAGGCAGTCTGGCTTCTCTCAGATGTGATTTGGAGACAAGAAGCAGCTCCTCACCCCTCAGGACCAATTCACAGGGCTC
760▶ ArgTrpGlnPheAsnArgGlnSerGlyPheLeuSerGlnMetTrpI leGlyAspLysLysGlnLeuLeuThrProLeuArgAspGlnPheThrArgAlaP
5201 CTCTGGACAATGACATTGGAGTGTCTGAGGCCACCAGGATTGACCCAAATGCTTGGGTGGAGAGGTGGAAGGCTGCTGGACACTACCAGGCTGAGGCTGC
793▶ roLeuAspAsnAspI leGlyValSerGluAlaThrArgI leAspProAsnAlaTrpValGluArgTrpLysAlaAlaGlyHisTyrGlnAlaGluAlaAl
5301 CCTGCTCCAGTGCACAGCAGACACCCTGGCTGATGCTGTCTGATCACCACAGCCATGCTTGGCAGCACAAGGCAAGACCCTGTTTCATCAGCAGAAAG
826▶ aLeuLeuGlnCysThrAlaAspThrLeuAlaAspAlaValLeuI leThrThrAlaHisAlaTrpGlnHisGlnGlyLysThrLeuPheI leSerArgLys
5401 ACCTACAGGATTGATGGCTCTGGACAGATGGCAATCACAGTGGATGTGGAGGTTGCCTCTGACACACCTCACCTGCAAGGATTGGCCTGAACTGTCAAC
860▶ ThrTyrArgI leAspGlySerGlyGlnMetAlal leThrValAspValGluValAlaSerAspThrProHisProAlaArgI leGlyLeuAsnCysGlnL
5501 TGGCACAGTGGCTGAGAGGGTGAAGTGGCTGGGCTTAGGCCCTCAGGAGAACTACCCTGACAGGCTGACAGCTGCCTGCTTTGACAGGTGGGACCTGCC
893▶ euAlaGlnValAlaGluArgValAsnTrpLeuGlyLeuGlyProGlnGluAsnTyrProAspArgLeuThrAlaAlaCysPheAspArgTrpAspLeuPr
5601 TCTGCTGACATGTACACCCTTATGTCTCCCTCTGAGAATGGCCTGAGGTGTGGCACCAGGAGCTGAACCTAGCTCACCAGTGGAGGAGAC
926▶ oLeuSerAspMetTyrThrProTyrValPheProSerGluAsnGlyLeuArgCysGlyThrArgGluLeuAsnTyrGlyProHisGlnTrpArgGlyAsp
5701 TTCCAGTTCACATCTCCAGTACTCTCAGCAACAGCTCATGGAACCTCTCACAGGCACCTGCCTCATGCAGAGGAGGAACCTGGCTGAACATTGATG
960▶ PheGlnPheAsnI leSerArgTyrSerGlnGlnGlnLeuMetGluThrSerHisArgHisLeuLeuHisAlaGluGluGlyThrTrpLeuAsnI leAspG
5801 GCTTCCACATGGGATGGAGGAGATGACTCTGGTCTCCTCTGTGTCTGCTGAGTTCAGTATCTGCTGGCAGGTACCCTATCAGCTGGTGTGGT
993▶ lyPheHisMetGlyI leGlyGlyAspSerTrpSerProSerValSerAlaGluPheGlnLeuSerAlaGlyArgTyrHisTyrGlnLeuValTrpCy

XbaI (5917)

5901 CCAGAAGTAAACCTAATCTAGAAGCTCGCTTTCTTGGCTGTCCAATTTCTATTAAGGTTCCCTTTGTTCCCTAAGTCCAACCTACTAACTGGGGATATTA
1026▶ sGlnLys•••
6001 TGAAGGGCCTTGAGCATCTGGATTCTGCCTAATAAAAAACATTTATTTTCATTGCAATGATGTATTTAAATTATTTCTGAATATTTTACTAAAAAGGAA
6101 TGTGGGAGGTCACTGCATTTAAAAACATAAAGAAATGAAGAGCTAGTTCAAACCTTGGGAAAATACACTATATCTTAAACTCCATGAAAGAAGGTGAGGCT
6201 GCAAAACAGCTAATGCACATTGGCAACAGCCCTGATGCCTATGCCTTATTCATCCCTCAGAAAAGGATTCAAGTAGAGGCTTGATTTGGAGGTTAAAGTT
PacI (6316)
6301 TTGCTATGCTGATTTTAAATTAATAAAACCCGCTTCGGCGGGTTTTTTTATGCATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCT
6401 TGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAAACCCGACAGGACTATAAAGATACCA
6501 GGGCTTTCCCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCGAGCCTGCGCTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTT
6601 TCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTCTGTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTAGCCGACCGCTGCGCCT
6701 TATCCGGTAACTATCGTCTTGGTCCAAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGG
6801 CGGTGTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAA
6901 AGAGTTGGTAGCTCTTGATCCGGCAAAACAAACCCGCTGGTAGCGGTGGTTTTTTTGTGGCAAGCAGCAGATTACCGCGCAAAAAAAGGATCTCAAG

7001 AAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGAGATTATCAAAAAGGATCTTCACCTA
7101 GATCCTTTTAAATTA AAAATGAAGTTTTAAATCAATCTAAAGTATATATGAGTAAACTTGGTCTGACAGTTACCAATGCTTAATCAGTGAGGCACCTATC
7201 TCAGCGATCTGTCTATTTTCGTTTCATCCATAGTTGCCTGACTCCCCGTCTGTAGATAACTACGATACGGGAGGGCTTACCATCTGGCCCCAGTGCTGCAA
7301 TGATACCGCGAGACCCACGCTCACCGGCTCCAGATTTATCAGCAATAAACAGCCAGCCGGAAGGGCCGAGCGCAGAAGTGGTCTGCAACTTTATCCGC
7401 CTCCATCCAGTCTATTAATTGTTGCCGGGAAGCTAGAGTAAGTAGTTCCGCCAGTTAATAGTTTGGCGAACGTTGTTGCCATTGCTACAGGCATCGTGGTG
7501 TCACGCTCGTCGTTTGGTATGGCTTCATTCAGCTCCGGTCCCAACGATCAAGGCGAGTTACATGATCCCCATGTTGTGCAAAAAAGCGTTAGCTCCT
7601 TCGGTCCTCCGATCGTTGTCAGAAGTAAGTTGGCCGAGTGTTATCACTCATGGTTATGGCAGCACTGCATAATTCTCTTACTGTCATGCCATCCGTAAG
7701 ATGCTTTTCTGTGACTGGTGAGTACTCAACCAAGTCACTCTGAGAATAGTGTATGCGGCAGCCGAGTTGCTCTTGCCCGGCTCAATACGGGATAATACC
7801 GCGCCACATAGCAGAACTTTAAAAGTGCTCATCATTGGAACGTTCTTCGGGGCGAAAACCTCAAGGATCTTACCGCTGTTGAGATCCAGTTCGATGT
7901 AACCCACTCGTGACCCAACTGATCTTCAGCATCTTTTACTTTTACCAGCGTTTCTGGGTGAGCAAAAACAGGAAGGCAAAAATGCCGAAAAAAGGGAAT
8001 AAGGGCGACACGAAATGTTGAATACTCATACTCTTCCTTTTTCAATATTATTGAAGCATTATCAGGGTATTGTCTCATGAGCGGATACATATTTGAA
8101 TGTATTTAGAAAAATAACAAATAGGGGTTCCGCGCACATTTCCCCGAAAAGTGCCACCTGACGTCTAAGAAACCATTATTATCATGACATTAACCTATA
8201 AAAATAGGCGTATCACGAGGCCCTTTCGTCTCGCGGTTTCGGTGATGACGGTGAAAACCTCTGACACATGCAGCTCCCGGAGACGGTCACAGCTTGTCT
8301 GTAAGCGGATGCCGGGAGCAGACAAGCCCGTCAAGGGCGGTCAGCGGGTGTGGCGGGTGTCCGGGCTGGCTTAACTATGCGGCATCAGAGCAGATTGTA
NdeI (8413)
8401 CTGAGAGTGCACCATATG