

pDUO-mTLR6/TLR2

A plasmid coexpressing the murine TLR6 and TLR2 genes

Catalog code: pduo-mtlr6tlr2

<https://www.invivogen.com/pduo-trl6-trl2>

For research use only

Version 20H25-MM

PRODUCT INFORMATION

Contents

- 20 µg of pDUO-mTLR6/TLR2 provided as 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

Toll-Like receptors (TLRs) play a critical role in early innate immunity to invading pathogens by sensing microorganisms. These evolutionary conserved receptors, homologues of the *Drosophila* Toll gene, recognize highly conserved structural motifs only expressed by microbial pathogens, called pathogen-associated microbial patterns (PAMPs). PAMPs include various bacterial cell wall components such as lipopolysaccharides (LPS), peptidoglycans and lipopeptides, as well as flagellin, bacterial DNA and viral double-stranded RNA. Stimulation of TLRs by PAMPs initiates a signaling cascade that involves a number of proteins, such as MyD88 and IRAK. This signaling cascade leads to the activation of the transcription factor NF-κB which induces the secretion of pro-inflammatory cytokines and effector cytokines that direct the adaptive immune response.

To date ten human and twelve murine TLRs have been characterized, TLR1 to TLR10 in humans, and TLR1 to TLR9, TLR11, TLR12 and TLR13 in mice, the homolog of TLR10 being a pseudogene. In many instances, TLRs require the presence of a co-receptor to initiate the signaling cascade. One example is TLR4 which interacts with MD2 and CD14 to induce NF-κB in response to LPS stimulation.

pDUO is an expression vector designed to co-express two TLRs or TLR-related genes known to interact with each other.

The genes cloned into pDUO comprise the coding sequence (without introns) from the ATG to the Stop codon.

PLASMID FEATURES

- **Murine TLR6** (2418 bp) / **Murine TLR2** (2352 bp)

TLR6 is expressed in spleen and BPL and, similarly to TLR1, acts as a co-receptor. Studies with dominant negative receptors have shown that TLR6 cooperates with TLR2 to recognize peptidoglycan and the yeast cell wall particle, zymosan¹. Furthermore, TLR6- and TLR2-deficient mice were reported to be hyporesponsive to mycoplasma macrophage-activating lipopeptide-2 kD (MALP-2), a diacylated lipoprotein, suggesting that TLR2 and TLR6 coordinate the response to this ligand. By contrast, TLR2 is able to recognize bacterial lipoproteins triacylated at the N-terminus cysteine residue². Thus TLR6 appears to discriminate between the N-terminal lipoylated structures of MALP-2 and lipopeptides derived from other bacteria.

- **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 protein³. 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. The enhancement varies from 2-fold in non-permissive cells to 20-fold in permissive cells. Furthermore, the SV40 enhancer is able to direct nuclear localization of plasmids⁴.

- **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 enhancer⁵.

- **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.*⁶

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

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

- **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 expression⁷.
- **EM7** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*.
- **Bsr (blasticidin resistance gene):** The *bsr* gene from *Bacillus cereus* encodes a deaminase that confers resistance to the antibiotic Blasticidin. In bacteria, *bsr* is expressed from the constitutive *E. coli* EM7 promoter. In mammalian cells, *bsr* is transcribed from the human FerH composite 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.

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 H₂O. 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.

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.

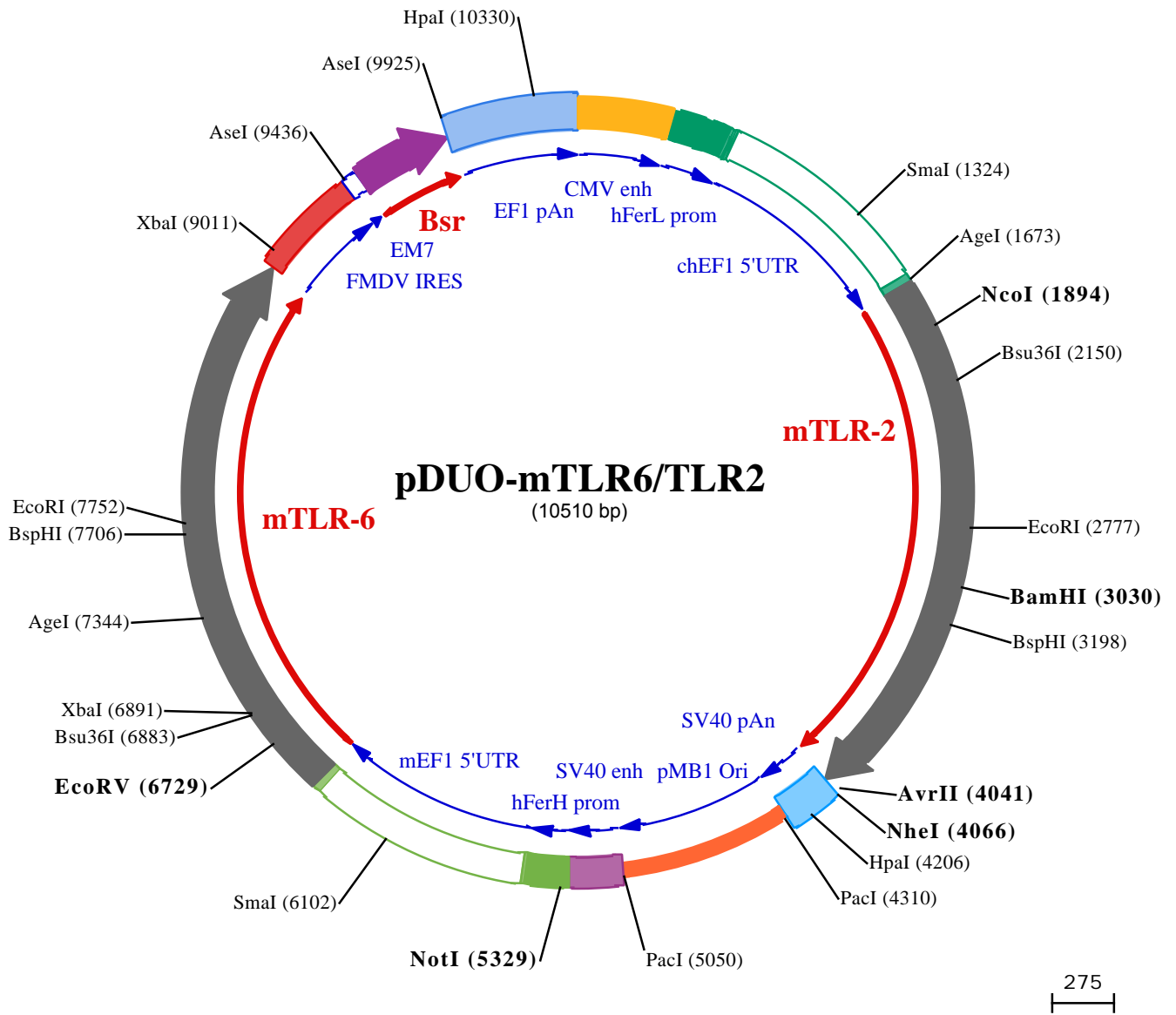
References

1. **Ozinsky A. et al., 2000.** The repertoire for pattern recognition of pathogens by the innate immune system is defined by cooperation between Toll-like receptors. *PNAS* 97(25):13766-71.
2. **Takeuchi O. et al., 2001.** Discrimination of bacterial lipoproteins by Toll-like receptor 6. *Int Immunol*, 13(7): 933-40.
3. **Eisenstein RS. & Munro HN. 1990.** Translational regulation of ferritin synthesis by iron. *Enzyme* 44(1-4):42-58.
4. **Dean DA. et al., 1999.** Sequence requirements for plasmid nuclear import. *Exp. Cell. Res.* 253:713-22.
5. **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.
6. **Carswell S. & Alwine JC. 1989.** Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. *Mol. Cell Biol.* 10: 4248-4258.
7. **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.

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 CCTGCAGCGTTACATAACTTACGGTAAATGGCCCGCTGGCTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAA
101 CGCCAATAGGGACTTTCCATTGACGTCAATGGGTGGAGTATTTACGGTAAACTGCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCC
201 TATTGACGTCAATGACGGTAAATGGCCCGCTGGCATTATGCCAGTACATGACCTTATGGGACTTTCTACTTGGCAGTACATCTACGTATTAGTCATC
301 GCTATTACCATGATGATCGGGTTTTGGCAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGGATTCCAAGTCTCCACCCATTGACGTCAATG
401 GGAGTTTGTCTTACTAGTCAGGCCCCAACCCCCCAAGCCCCATTTCACAACAGCTGGCGCTACAGGCGCTGACTTCCCCCTGCTTTGGGGCGGG
501 GGGCTGAGACTCTATGTGCTCCGGATTGGTCAGGCAGGCGCTTCGGCCCGCTCCTGCCACGCGAGATTGCGCGTAGCCTCCCGAGCGCCCTGCC
601 TCCGAGGCGCGGCACCATAAAGAAGCCGCCCTAGCCACGTCCTCCCGCAGTTCGGCGGTCCCGGGTCTGTCTCAAGCTTGGCGCAGAACACAGg
701 taagtgccgtgtgtggtccccgggctggcctctttacgggttatggccttgcgtgcctgaattacttccatgccctggctgcagtacgtgatc
801 ttgatcccagacttccgggttgggaagtgggtgggagagttcgaggccttgcgcttaaggagcccttgcctcgtgcttgagttgaggcctggcttgggagc
901 ctggggcccgcgctgctaactcgggtggcaccttcgcgctgtctcgctgcttgcctaaagtctctagccatttaaattttgat aaccagctgcgacg
1001 cttttttctggcgagatagctcttgaatcggggccaggatctgcacactggtatctcggttttggggccgaggcgaggggcccgtgcgtccc
1101 agcgcacatgctcggcgaggcggggcctgcgagcgcggccaccgagaatcgacgggggtagctctcaactggccgctgctctggtgcctggcctcgc
1201 gccccgctgtatgccccgcccctgggcccgaaggctggcccggctcgccaccagttgcgtgagcggaaagatggccgcttcccggcctgctgcaggagc

SmaI (1324)

1301 tcaaaatggaggacgcccgggagagcggggcgggtgagtcacccacacaaaggaaaaggccttctctcatccgctgcttcatgtgactcca
1401 cggagtagccggcgccgtccaggcacctcgattagttgtcgagctttggagtagctcgtctttaggttgggggaggggttttatgcatggagttcc
1501 ccacactgagtggtggagactgaagagttaggccagcttggcacttgatgtaattctccttggaaattgccttttgagttggatcttgcctcattc

AgeI (1673)

1601 tcaagcctcagacagtggttcaaagttttttcttccatttcagGTGTCGTGAAAACCTACCCCTAAAAGCCCGTAGGAGGGCCAGCATGCTACGAGC
1701 TCTTTGGCTCTTCTGGATCTTGGTGGCCATAACAGTCTCTTCAGCAAACGCTGTTCTGCTCAGGAGTCTCTGTCTATGTGATGCTTCTGGGGTGTGTGAT
4 aLeuTrpLeuPheTrpI leLeuValAlaI leThrValLeuPheSerLysArgCysSerAlaGlnGluSerLeuSerCysAspAlaSerGlyValCysAsp

NeoI (1894)

1801 GGCCGCTCCAGGTCTTTACCTCTATTCCCTCCGACTCACAGCAGCCATGAAAAGCCTTGACTGTCTTTCAACAAGATCACCTACATTGGCCATGGTG
38 GlyArgSerArgSerPheThrSerI leProSerGlyLeuThrAlaAlaMetLysSerLeuAspLeuSerPheAsnLysI leThrTyrI leGlyHisGlyA
1901 ACCTCGAGCGTGTGCGAACCTCCAGTTCTGATTTGAAGTCCAGCAGAATCAATACAATAGAGGGAGACGCTTTTATTCTCTGGGCAGTCTTGAACA
71 spLeuArgAlaCysAlaAsnLeuGlnValLeuI leLeuLysSerSerArgI leAsnThrI leGluGlyAspAlaPheTyrSerLeuGlySerLeuGluHi
2001 TTTGGATTTGCTGATAATCACCTATCTAGTTTATCTTCTCTGGTTCGGGCCCTTCTCTTTGAAATACTTAAACTTAATGGGAAATCCTTACCAG
104 sLeuAspLeuSerAspAsnHisLeuSerSerLeuSerSerTrpPheGlyProLeuSerSerLeuLysTyrLeuAsnLeuMetGlyAsnProTyrGln

Bsu36I (2150)

2101 AACTGGGGTAAACATCGCTTTTCCCAATCTCACAAATTTACAAAACCTCAGGATAGGAAATGTAGAGACTTTCAGTGAGATAAGGAGAATAGATTTTG
138 ThrLeuGlyValThrSerLeuPheProAsnLeuThrAsnLeuGlnThrLeuArgI leGlyAsnValGluThrPheSerGluI leArgArgI leAspPheA
2201 CTGGGCTGACTTCTCAATGAACCTTGAATTAAGGCATTAAGTCTCCGGAATTATCAGTCCCAAAGTCTAAAGTCGATCCGCGACATCCATCAGCTGAC
171 laGlyLeuThrSerLeuAsnGluLeuGluI leLysAlaLeuSerLeuArgAsnTyrGlnSerGlnSerLeuLysSerI leArgAspI leHisHisLeuTh
2301 TCTTCACTTAAGCGAGTCTGCTTTCCCTGCTGGAGATTTTTCAGATATTTCTGAGTTCTGTGAGATATTTAGAACTAAGAGATACTAAGTGGCCAGGTT
204 rLeuHisLeuSerGluSerAlaPheLeuLeuGluI lePheAlaAspI leLeuSerSerValArgTyrLeuGluLeuArgAspThrAsnLeuAlaArgPhe
2401 CAGTTTCCACACTGCCGTAGATGAAGTCAGCTCAGCATCCGATGAAGAAGTCCGAGGCTCGGTTCTCACTGATGAAAGCTTAAACGAGCTCCTGA
238 GlnPheSerProLeuProValAspGluValSerSerProMetLysLysLeuAlaPheArgGlySerValLeuThrAspGluSerPheAsnGluLeuLeuL
2501 AGCTGTTGCGTTACATCTTGAAGTCTCGGAGGTAGAGTTCGACGACTGTACCTCAATGGGCTCGGCGATTTCACCCCTCGGAGTCAGACGTAGTGAG
271 ysLeuLeuArgTyrI leLeuGluLeuSerGluValGluPheAspAspCysThrLeuAsnGlyLeuGlyAspPheAsnProSerGluSerAspValValSe
2601 CGAGCTGGTAAAGTAGAAACAGTCACTATCCGAGGTTGCATATCCCCAGTCTATTTGTTTTATGACCTGAGTACTGTCTATCCCTCCTGGAGAAG
304 rGluLeuGlyLysValGluThrValThrI leArgArgLeuHisI leProGlnPheTyrLeuPheTyrAspLeuSerThrValTyrSerLeuLeuGluLys

EcoRI (2777)

2701 GTGAAGCAATCAGAGTACAGAACAGCAAGTCTTCTGCTGCTGCTTCTCCAGCATTTAAAATCATTAGAAATCTTAGACCTCAGCGAAAATC
338 ValLysArgI leThrValGluAsnSerLysValPheLeuValProCysSerPheSerGlnHisLeuLysSerLeuGluPheLeuAspLeuSerGluAsnL
2801 TGATGGTTGAAGAATATTGAAGAACCTCAGCTGTAAAGGAGCCTGGCCTTCTCTACAAACCTAGTTTGTAGCCAGAATCATTGAGATCAATGCAAAA
371 euMetValGluGluTyrLeuLysAsnSerAlaCysLysGlyAlaTrpProSerLeuGlnThrLeuValLeuSerGlnAsnHisLeuArgSerMetGlnLy
2901 AACAGGAGAGATTTGCTGACTCTGAAAAACCTGACCTCCCTTGACATCAGCAGGAACACTTTTCATCCGATGCCGACAGCTGTGAGTGGCCAGAAAAG
404 sThrGlyGluI leLeuLeuThrLeuLysAsnLeuThrSerLeuAspI leSerArgAsnThrPheHisProMetProAspSerCysGlnTrpProGluLys

BamHI (3030)

3001 ATGCGCTTCTGAATTTGTCCAGTACAGGGATCCGGGTGGTAAAAACGTCATTCTCAGACGCTGGAGGTGTTGGATGTTAGTAAACAATCTTGACT
438 MetArgPheLeuAsnLeuSerSerThrGlyI leArgValValLysThrCysI leProGlnThrLeuGluValLeuAspValSerAsnAsnLeuAspS

BspHI (3198)

3101 CATTTTCTTGTCTTGGCTCGGCTGCAAGAGCTCTATATTTCCAGAAATAAGCTGAAAACACTCCAGATGCTTCGTTGTTCCCTGTGTTGCTGGTCAT
471 erPheSerLeuLeuProArgLeuGlnGluLeuTyrI leSerArgAsnLysLeuLysThrLeuProAspAlaSerLeuPheProValLeuLeuValMe
3201 GAAAAATCAGAGAGAATGCAGTAAAGTACTTTCTTAAAGACCAACTTGGTTCTTTTCCAAAACCTGGAGACTCTGGAAGCAGCGACAACCACTTTGTTTGC
504 tLysI leArgGluAsnAlaValSerThrPheSerLysAspGlnLeuGlySerPheProLysLeuGluThrLeuGluAlaGlyAspAsnHisPheValCys
3301 TCCTGCGAACTCCTATCCTTTACTATGGAGACGCCAGCTCTGGCTCAAATCCTGGTTGACTGGCCAGACAGTACCTGTGTGACTCTCCGCTCGCCTGC
538 SerCysGluLeuLeuSerPheThrMetGluThrProAlaGlnI leLeuValAspTrpProAspSerTyrLeuCysAspSerProArgLeuH
3401 ACGCCACAGGCTTACAGATGCCCGCCCTCGTCTTGAATGTACCAAGCTGCAGCTGGTGTCTGGAGTCTGCTGCGCCTTCTCTGTTGATCTGTCT
571 isGlyHisArgLeuGlnAspAlaArgProSerValLeuGluCysHisGlnAlaAlaLeuValSerGlyValCysCysAlaLeuLeuLeuLeuI leLeuLe

3501 CGTAGGTGCCCTGTGCCACCATTCCACGGACTGTGGTACCTGAGAATGATGTGGGCGTGGCTCCAGGCCAAGAGGAAGCCCAAGAAAGCTCCCTGCAGG
604 uValGlyAlaLeuCysHisHisPheHisGlyLeuTrpTyrLeuArgMetMetTrpAlaTrpLeuGlnAlaLysArgLysProLysLysAlaProCysArg
3601 GACGTTTTGCTATGATGCCTTTGTTTCACAGTGCAGGATTCCCATTTGGGTGGAGAACCCTCATGGTCCAGCAGCTGGAGAACCCTGACCCGCCCTTTA
638 AspValCysTyrAspAlaPheValSerTyrSerGluGlnAspSerHisTrpValGluAsnLeuMetValGlnGlnLeuGluAsnSerAspProProPheL
3701 AGCTGTGTCTCCACAAGCGGGACTTCGTTCCGGGCAAAATGGATCATTGACAACATCATCGATTCCATCGAAAAGAGCCACAAAACCTGTGTTCGTGCTTTC
671 ysLeuCysLeuHisLysArgAspPheValProGlyLysTrpIleIleAspAsnIleIleAspSerIleGluLysSerHisLysThrValPheValLeuSe
3801 TGAGAACCCTCGTACGGAGCGAGTGGTCAAAGTACGAACCTGGACTTCCCACTTCAGGCTCTTGACGAGAACAACGACGGCCATCCTGTTTTGCTG
704 rGluAsnPheValArgSerGluTrpCysLysTyrGluLeuAspPheSerHisPheArgLeuPheAspGluAsnAsnAspAlaAlaIleLeuValLeuLeu
3901 GAGCCATTGAGAGGAAAGCCATTCCCAGCGCTTCTGCAAACTGCGCAAGATAATGAACACCAAGACCTACCTGGAGTGGCCCTGGATGAAGGCCAGC
738 GluProlIleGluArgLysAlaIleProGlnArgPheCysLysLeuArgLysIleMetAsnThrLysThrTyrLeuGluTrpProLeuAspGluGlyGlnG

AvrII (4041)

NheI (4066)

4001 AGGAAGTGTTTTGGGTAATCTGAGAACTGCAATAAAGTCCTAGTTCCTCCACCCAGTTCCTGAGCTAGCTGGCCAGACATGATAAGATACATTGATGAG
771 InGluValPheTrpValAsnLeuArgThrAlaIleLysSer...
4101 TTTGGACAAACCACTAGAATGCAGTGAAAAAATGCTTTATTGTGAAATTGTGATGCTATTGCTTTATTGTAAACCATTATAAGCTGCAATAAAC

HpaI (4206)

4201 AAGTTAAACAACAACAAATTGCATTCTTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTTAAAGCAAGTAAACCTCTACAAATGTGTTATGGA

PacI (4310)

4301 AATGTTAATTAACATAGCCATGACCAAAATCCCTTAACGTGAGTTTTTCGTTCCACTGAGCGTCAGACCCCGTAGAAAAGATCAAAGGATCTTCTTGAGATC

4401 CTTTTTTTTCTGCGCGTAATCTGCTGCTTGAACAAAAAACCACCGCTACCAGCGGTGTTTGTGTTGCCGATCAAGAGCTACCAACTCTTTTTCCGAA

4501 GGTAACGGCTTCAGCAGAGCGCAGATACCAAACTGTTCTTCTAGTGTAGCCGATGTTAGGCCACCACTCAAGAAGCTGTAGCACCAGCTACATAC

4601 CTCGCTCTGCTAATCTGTTACCAGTGGCTGCTGCCAGTGGCGATAAGTCTGTCTTACCAGGTTGGACTCAAGACGATAGTTACCAGGATAAGGCGCAGC

4701 GGTCCGGCTGAACGGGGGTTCTGTGCACACAGCCAGCTGGAGCGAACGACCTACACCGAAGTGGAGATACCTACAGCGTGGACTATGAGAAAGCGCCAC

4801 GCTTCCCAAGGGAGAAAGGCGGACAGGTATCCGGTAAGCGGCAGGGTCCGAACAGGAGAGCGCAGGAGGAGCTTCCAGGGGAAACGCTGGTATCTT

4901 TATAGTCTGTGGGTTTCGCCACCTCTGACTTGAGCGTGCATTTTTGTGATGCTCGTCAGGGGGCGGAGCCTATGGAAAAACGCCAGCAACCGCGCCT

PacI (5050)

5001 TTTTACGGTTCCTGGCCTTTTGTGCGCTTTTGTCTCACATGTTCTTAATTAACCTGCAGGGCCTGAAATAACCTCTGAAAGAGGAACTTGGTTAGGTACC

5101 TTCTGAGGCTGAAAGAACCAGCTGTGGAATGTGTGTCAGTTAGGGTGTGGAAAGTCCCAGGCTCCCAGCAGGCAGAAATGCAAAGCATGCATCTCA

5201 ATTAGTCAGCAACCAGGTGTGAAAGTCCCAGGCTCCCAGCAGGCAGAAATGCAAAGCATGCATCTCAATTAGTCAGCAACCATAGTCCCACTAGT

NotI (5329)

5301 TCCGCCAGAGCGCGAGGGCCTCCAGCGCGCCCTCCCCACAGCAGGGGGGGGTCCCGCGCCACCGGAAGGAGCGGGCTCGGGCGGGCGGGCGC

5401 TGATTGGCCGGGGCGGGCTGACGCCGACCGGCTATAAGAGACCACAAGCAGCCCGCAGGGCCAGACGTTCTTCGCCGAAGCTTGCCCTCAGAACGCAG

5501 gtgaggggcggggtgtggcttccgcgggccgagctggaggtcctgctccgagcgggccccgctgtcgctcggggggattagctgcgagcattc

5601 ccgcttcgagttcgggcgggcggggagcagagtgagggcctagcggcaaccccgtagcctcgctcggttcgagccttagggcctagcgtggtgtccg

5701 cgccgcccgcgctgctactccggccgactctggtctttttttttttgtgtgtgctgctgcttcgattgccgttcagcaataggggctaaca

5801 aagggagggtgccccgcttgcctcgcccggagcccggagaggtcatggttggggaggaatggagggacaggagtgccggctggggcccccgccctcgga

5901 gcacatgcccagccacactggatggggcgaggcctggggttttcccgaagcaaccaggctggggttagcgtgcccaggccatgtggccccagcaccg

SmaI (6102)

6001 gcacgatctggcttggcggcgccgcttgcctgcctcctaactagggtgagccatcccgtccggcaccagttgcgtgctggaagatggccgctcc

6101 cgggccctgttgcaaggagctcaaaatggaggacgcccagcagccggaggcggggtgagtcacccacaaaaggaagggcctggtccctcaccg

6201 gctgctgcttccctgtgaccccggtcctatcgcccgcaatagtcacctcgggcttttgagcacggctagtcgcccggggggaggggatgtaatggcgt

6301 tggagtttggttcacatttgggtgggtggagactagtcaggccagcctggcgtggaagtcattttggaaattgtcccttgagtttgagcggagctaat

6401 tctcgggcttcttagcgggttcaaaggtatcttttaacccttttttagGTGTTGTGAAAACCACCGCTAATTCAAAGCAATCATGGTAAAGTCCCTCTGG

MetValLysSerLeuTrp

6501 GATAGCCTCTGCAACATGAGCCAAGACAGAAAACCCATCGTGGGAGTTTCCACTTTGTTTGCGCCCTGGCCTTAATAGTCGGAAGCATGACCCGTTCT

7 AspSerLeuCysAsnMetSerGlnAspArgLysProIleValGlySerPheHisPheValCysAlaLeuAlaLeuIleValGlySerMetThrProPheS

6601 CTAATGAACCTGAGTCTATGTTAGACTATTCACACAGGAACCTTACTCATGTCCCCAAAGACCTGCCACCAAGAACAAAAGCCCTGAGTCTGTCTCAAAA

40 erAsnGluLeuGluSerMetValAspTyrSerAsnArgAsnLeuThrHisValProLysAspLeuProProArgThrLysAlaLeuSerLeuSerGlnAs

EcoRV (6729)

6701 CTCTATATCTGAGCTTCGGATGCCTGATATCAGCTTCTGTCTCAGAGCTGAGAGTCTGAGACTCTCCCAACACAGGATACGGAGCCTTGATTTCCATGTA

73 nSerIleSerGluLeuArgMetProAspIleSerPheLeuSerGluLeuArgValLeuArgLeuSerHisAsnArgIleArgSerLeuAspPheHisVal

XbaI (6891)

Bsu36I (6883)

6801 TTCTGTTCATCAGACTTAGAATACCTGGATGTCTCACAAATCGGTTGCAAAACATCTCTGTGTCGCCCTATGGCGAGCCTGAGGCATCTAGACCTCT

107 PheLeuPheAsnGlnAspLeuGluTyrLeuAspValSerHisAsnArgLeuGlnAsnIleSerCysCysProMetAlaSerLeuArgHisLeuAspLeuS

6901 CATTCAATGACTTGTGACTGCCTGTGTAAGGAATTTGGCAACCTGACGAAGCTGACTTCTCGGATTAAGTGTGCAAGTTCGCAACACTGGA

140 erPheAsnAspPheAspValLeuProValCysLysGluPheGlyAsnLeuThrLysLeuThrPheLeuGlyLeuSerAlaAlaLysPheArgGlnLeuAs

7001 TCTGCTCCAGTTGCTCACTTGCATCTAAGCTGCATTCTTCTGACTTAGTGAGTCATCATATAAAAGCGGGGAAACAGAAAGTCTTCAGATTCCCAAT
173▶ pLeuLeuProValAlaHisLeuHisLeuSerCysI leLeuLeuAspLeuValSerHisHisI leLysGlyGlyGluThrGluSerLeuGlnI leProAsn
7101 ACCACCGTTCTCCATTTGGTCTTTCATCCAAAATAGCTTGTCTCTGTTCAAGTGAACATGTCTGTAAACGCTTTAGGACATTTACAACCTGAGTAATATTA
207▶ ThrThrValLeuHisLeuValPheHisProAsnSerLeuPheSerValGlnValAsnMetSerValAsnAlaLeuGlyHisLeuGlnLeuSerAsnI leL
7201 AATTGAATGATGAAAAGTCAAGGTTAATGACATTTTATCAGAACTCACAGAGGTCACACCTTATTGAATGTGACCCCTCAGCACATAGAAAACAAC
240▶ ysLeuAsnAspGluAsnCysGlnArgLeuMetThrPheLeuSerGluLeuThrArgGlyProThrLeuLeuAsnValThrLeuGlnHisI leGluThrTh

AgeI (7344)

7301 CTGGAAGTGTCTGGTAAACTTTTCCAATTCTTTTGGCCCCGACCGGTGGAGTACCTCAATATTTACAACCTTAAACGATAACTGAGAGAATCGACAGGGAA
273▶ rTrpLysCysSerValLysLeuPheGlnPhePheTrpProArgProValGluTyrLeuAsnI leTyrAsnLeuThrI leThrGluArgI leAspArgGlu
7401 GAATTTACTTACTCGGAGACAGCACTGAAGTCACTGATGATAGAGCAGCTCAAAAACCAAGTGTCTCTTTTCAAAGGAGGCGCTATACCTGGTGTGTTG
307▶ GluPheThrTyrSerGluThrAlaLeuLysSerLeuMetI leGluHisValLysAsnGlnValPheLeuPheSerLysGluAlaLeuTyrSerValPheA
7501 CTGAGATGAACATCAAGATGCTCTCTATCTCAGACACCCCTTTTATCCACATGGTGTGCCGCCATCCCCAAGCTCATTTACATTTCTGAACCTTACCCA
340▶ laGluMetAsnI leLysMetLeuSerI leSerAspThrProPheI leHisMetValCysProProSerProSerSerPheThrPheLeuAsnPheThrGl
7601 GAATGTTTTTACTGACAGTGTGTTTCAAGGCTGTCCACCTTAAAGAGATTGCAGACACTTATCTTACAAGGAATGGTTTGAAGAACTTTTTTAAAGTA
373▶ nAsnValPheThrAspSerValPheGlnGlyCysSerThrLeuLysArgLeuGlnThrLeuI leLeuGlnArgAsnGlyLeuLysAsnPhePheLysVal

BspHI (7706)

EcoRI (7752)

7701 GCTCTCATGACTAAGAATATGCTCTCTCTGAAACTTTGGATGTTAGTTTGAATCTTTGAACTCTCATGCATATGACAGGACATGCGCCTGGGCTGAGA
407▶ AlaLeuMetThrLysAsnMetSerSerLeuGluThrLeuAspValSerLeuAsnSerLeuAsnSerHisAlaTyrAspArgThrCysAlaTrpAlaGluS
7801 GCATATTGGTGTGTAATTTGCTTCCGAATATGCTTACAGGCTCTGTCTTCCAGATGCTTACCTCCCAAGTCAAGGTCCTTGACCTTCAACAACAGGAT
440▶ erI leLeuValLeuAsnLeuSerSerAsnMetLeuThrGlyValPheArgCysLeuProLysValLysValLeuHisGlnAsnArgI I
7901 AATGACCTCCCTAAAGATGTCAACCCAGCTGCAGGCTTTGCAAGAACTCAATGTAGCATCCAACCTCTTAACTGACCTTCCGGGTGGGGCCTTACGC
473▶ eMetSerI leProLysAspValThrHisLeuGlnAlaLeuGlnGluLeuAsnValAlaSerAsnSerLeuThrAspLeuProGlyCysGlyAlaPheSer
8001 AGCCTTTCTGTGCTGCTCATCGACCATAACTCAGTTTCCCATCCCTCTGAGGATTTCTTCCAGAGCTGTGAGAAATATTAGATCCCTAACAGCGGGAACA
507▶ SerLeuSerValLeuVal I leAspHisAsnSerValSerHisProSerGluAspPhePheGlnSerCysGlnAsnI leArgSerLeuThrAlaGlyAsnA
8101 ACCCATTTCCAATGCACATGTGAGCTGAGGGACTTTGTCAAGAACATAGGCTGGGTAGCAAGAGAAGTGGTGGAGGCGTGGCCTGACTCTTGAACGTTGTA
540▶ snProPheGlnCysThrCysGluLeuArgAspPheValLysAsnI leGlyTrpValAlaArgGluValValGluGlyTrpProAspSerTyrArgCysAs
8201 CTACCAGAAAGCTTAAGGGAAGTGCAGTGCAGGGACTTCCACATGCTCCACTGCTGTGACTGTTCTGTGACTGTACCATCGGGGCCACTATG
573▶ pTyrProGluSerSerLysGlyThrAlaLeuArgAspPheHisMetSerProLeuSerCysAspThrValLeuLeuThrValThrI leGlyAlaThrMet
8301 CTGGTGTGGCTGCTACTGGGGCTTTCCTCTGTCTACTTTGACCTGCCCTGGTATGTGAGGATGTGTGTGCTGAGTGGACACAGACCAGGCACAGGGCCA
607▶ LeuValLeuAlaValThrGlyAlaPheLeuCysLeuTyrPheAspLeuProTrpTyrValArgMetLeuCysGlnTrpThrGlnThrArgHisArgAlaA
8401 GGCACATCCCCTTAGAGGAACTCCAGAGAACTCCAGTCCATGCTTTTGTCTCATAACAGTGCAGTATGTTCTGCCTGGGTGAGAAACGAATTACTACC
640▶ rgHisI leProLeuGluGluLeuGlnArgAsnLeuGlnPheHisAlaPheValSerTyrSerGluHisAspSerAlaTrpValLysAsnGluLeuLeuPr
8501 CAACCTAGAGAAAGATGACATCCGGGTTTGCCTCCATGAGAGGAACCTTTGTCCCTGGCAAGAGCATTGTGGAGAATCATCAATTTTATTGAGAAGAGT
673▶ oAsnLeuGluLysAspAspI leArgValCysLeuHisGluArgAsnPheValProGlyLysSerI leValGluAsnI leI leAsnPheI leGluLysSer
8601 TACAAGGCCATCTTTGTGCTGTCTCCCACTTCATCCAGAGTGAGTGGTCCATTATGAACTCTATTTTGGCCATCATAATCTTCCATGAAGGCTCTG
707▶ TyrLysAlaI lePheValLeuSerProHisPheI leGlnSerGluTrpCysHisTyrGluLeuTyrPheAlaHisHisAsnLeuPheHisGluGlySerA
8701 ATAACCTAATCTCATCTTGTGGAACCCATTTACAGAACAACATTTCCAGTAGATACCAAGCTGCGGGCTCTCATGGCACAGCGGACTTACTTGGGA
740▶ spAsnLeuI leLeuI leLeuLeuGluProI leLeuGlnAsnAsnI leProSerArgTyrHisLysLeuArgAlaLeuMetAlaGlnArgThrTyrLeuGI
8801 ATGGCTACTGAGAAGGGCAAACGTTGGGCTGTTTGGGCCAACCTTAGAGCTTCATTTATTATGAAGTTAGCCTTAGTCAATGAGGATGATGTGAAAAC
773▶ uTrpProThrGluLysGlyLysArgGlyLeuPheTrpAlaAsnLeuArgAlaSerPheI leMetLysLeuAlaLeuValAsnGluAspAspValLysThr
8901 TGAACCTTGGGTTTCTAACTTAATAAATCTGCAACCTGGGCTCTGTAGGAGCAGGTTTCCCAATGACACAAAACCTGCAACTTGAACCTCCGCTGGT
807▶ ...

XbaI (901)

9001 CTTTCCAGGCTAGAGGGTAACACTTTGTACTGCGTTTGGCTCCACGCTCGATCCACTGGCGAGTGTAGTAAACAGCACTGTTGCTTCGTAGCGGAGCA
9101 TGACGGCCGTGGGAACTCCTCCTTGTAACAAGGACCCACGGGGCCAAAAGCCACGCCCACAGGGCCCGTCTGTGTGCAACCCAGCAGCGGGACTTT
9201 ACTGCGAAACCCACTTTAAAGTGACATTGAAACTGGTACCCACACTGGTGACAGGCTAAGGATGCCCTTACAGTACCCCGAGGTAACACGCGCACTC
9301 GGGATCTGAGAAGGGGACTGGGGCTTCTATAAAAGCGCTCGGTTTAAAAAGCTTCTATGCCTGAATAGGTGACCGGAGGTGGCACCTTTCTTTGCAAT

AseI (9436)

9401 TACTGACCTATGAATACACTGACTGTTTGACAATTAATCATCGGCATAGTATATCGGCATAGTATAATACGACTCACTATAGGAGGGCCACCATGAAGA
9501 CCTTCAACATCTCTCAGCAGGATCTGGAGCTGGTGGAGGTCGCCACTGAGAAGATCACCATGCTCTATGAGGACAACAAGCACCATGTGGGGCGGCCAT
3▶ hrPheAsnI leSerGlnGlnAspLeuGluLeuValGluValAlaThrGluLysI leThrMetLeuTyrGluAspAsnLysHisHisValGlyAlaAlaI
9601 CAGGACCAAGACTGGGAGATCATCTCTGTGTCCACATTGAGGCTACATTGGCAGGGTCACTGTCTGTGCTGAAGCCATTGCCATTGGGTCTGTGTG
36▶ eArgThrLysThrGlyGluI leI leSerAlaValHisI leGluAlaTyrI leGlyArgValThrValCysAlaGluAlaI leAlaI leGlySerAlaVal
9701 AGCAACGGGCAGAAAGACTTTGACACCATTTGTGGCTGTGAGCACCCTACTCTGATGAGGTGGACAGATCCATCAGGGTGGTCAGCCCCTGTGGCATGT
70▶ SerAsnGlyGlnLysAspPheAspThrI leValAlaValArgHisProTyrSerAspGluValAspArgSerI leArgValValSerProCysGlyMetC
9801 GCAGAGAGCTCATCTGACTATGCTCCTGACTGCTTGTGTCTATTGAGATGAATGGCAAGCTGGTCAAAACCCACCATTTAGGAACTCATCCCCCTCAA
103▶ ysArgGluLeuI leSerAspTyrAlaProAspCysPheValLeuI leGluMetAsnGlyLysLeuValLysThrThrI leGluGluLeuI leProLeuLy

AseI (9925)

9901 GTACACCGAAGAACTAAACCTGAATTAATTCGTAGGATTATCCCTAATACCTGCCACCCCACTTAACTCAGTGGTGAAGAACGGTCTCAGAAGTGT
136▶ sTyrThrArgAsn...
10001 GTTTCAATTTGGCCATTTAAGTTTAGTAGTAAAAGACTGGTTAATGATAACAATGCATCGTAAAACCTTCAAGAAAGGAGAATGTTTTGTGGACCACT
10101 TTGGTTTTCTTTTTGCGTGTGGCAGTTTAAAGTTATTAGTTTTTAAAAATCAGTACTTTTTAATGGAAAACACTTGACCAAAAAATTTGTACAGAATTTT
10201 GAGACCCATTAATAAAGTTAAATGAGAAACCTGTGTGTTCTTTGGTCAACACCCAGACATTTAGTGAAAGACATCTAATTTCTGGTTTTACGAATCTGG

HpaI (10330)

10301 AAACCTCTTGAAAATGTAATCTTGTAGTTAACACTTCTGGGTGGAGAATAGGTTGTTTTCCCCCACATAATTGGAAGGGGAAGGAATATCATTAAAG
10401 CTATGGGAGGTTTCTTTGATTACAACACTGGAGAGAAATGCAGCATGTTGCTGATTGCCTGTCACTAAAACAGGCCAAAAACTGAGTCTTGGGTTGCA
10501 TAGAAAGCTG