

# pDUO-mTLR1/TLR2

A plasmid coexpressing the murine TLR1 and TLR2 genes

Catalog code: pduo-mtlr1tlr2

<https://www.invivogen.com/pduo-trl1-trl2>

For research use only

Version 20H25-MM

## PRODUCT INFORMATION

### Contents

- 20 µg of pDUO-mTLR1/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 TLR1 (2388 bp) / Murine TLR2 (2355 bp)**

TLR1, the first member of the Toll-Like Receptor family was identified by the presence of a domain homology found in both *Drosophila* Toll and human IL-1 receptors. So far, no direct ligands for TLR1 have been identified however it seems to act as a coreceptor. TLR1 has been shown to associate with TLR2 in response to triacylated lipopeptides<sup>1</sup>, but not diacylated lipopeptides<sup>2</sup>. These observations suggest that TLR1 is able to discriminate among lipoproteins by recognizing the lipid configuration.

- **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 protein<sup>3</sup>. 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 plasmids<sup>4</sup>.

- **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 enhancer<sup>5</sup>.

- **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.*<sup>6</sup>

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

- **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 expression<sup>7</sup>.

## 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](mailto:info@invivogen.com)



- 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 H<sub>2</sub>O. 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

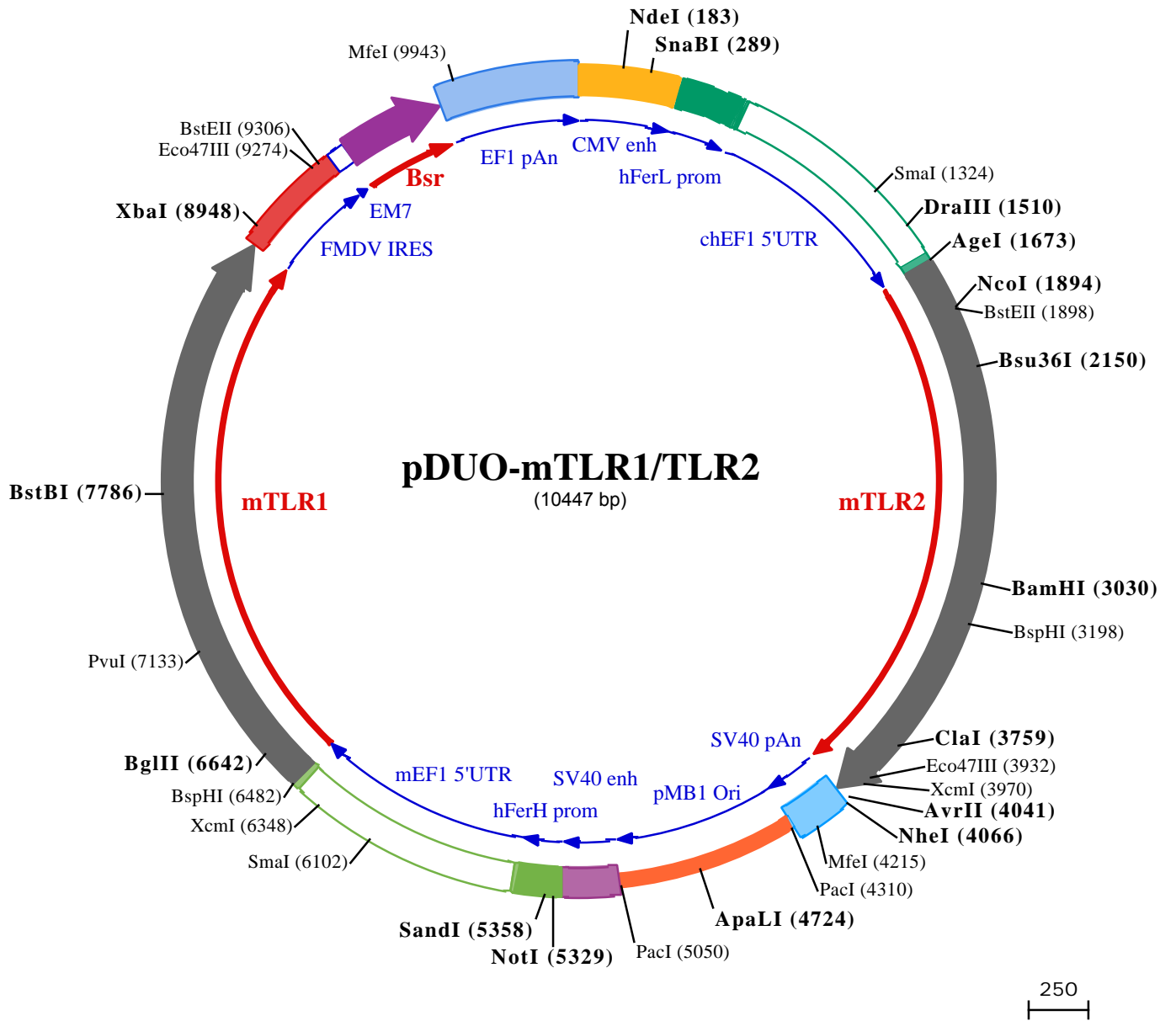
1. **Takeuchi O. et al., 2002.** Cutting edge: Role of Toll-like receptor 1 in mediating immune response to microbial lipoproteins. *J Immunol*, 169(1):10-14.
2. **Takeuchi O. et al., 2001.** Discrimination of bacterial lipoproteins by Toll-like receptor 6. *Int Immunol*, 13(7): 933-40.
3. **Eisenstein 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](mailto:info@invivogen.com)





1 CCTGCAGCGCTTACATAACTTACGGTAAATGGCCCGCTGGCTGACCGCCCAACGACCCCGCCATTGACGTCATAATGACGTATGTTCCCATAGTAA

NdeI (183)

101 CGCCAAATAGGGACTTTCATTGACGTCAATGGGTGGAGTATTTACGGTAAACTGCCCACTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCC

SnaBI (289)

201 TATTGACGTCAATGACGGTAAATGGCCCGCTGGCATTATGCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATC

301 GCTATTACCATGATGATGCGGTTTTGGCAGTACATCAATGGGCGTGATAGCGGTTTACTACGGGGATTCCAAGTCTCCACCCATTGACGTCAATG

401 GGAGTTTGTTTGACTAGTCAGGGCCCAACCCCCCAAGCCCCATTTACAACACGCTGGCGCTACAGGCGGTGACTTCCCCTTGCTTTGGGGCGGG

501 GGGCTGAGACTCCTATGTGCTCCGATTGGTCAGGCAGGCGCTTCGGCCCCGCTCTGCCACCGAGATTGGCCGTAGGCCCTCCCCGAGCCCTGCC

601 TCCGAGGGCCGGCCACCATAAAAAGAGCCGCCCTAGCCACGTCCTCCGAGTTCGGCGGTCCCAGGGTCTGTCTAAAGCTTGCCGCCAGAACACAGG

701 taagtgccgtgtgtggttcccggggcctggcctctttacgggttatggccttgcgtgccttgaattacttccatgccctggctgcagtacgtgatctc

801 ttgatcccagccttcggggttgaagtgggtgggagagttcgaggccttgcgcttaaggagccccttgcctcgtgcttgagttgaggcctggcttgggcg

901 ctggggcccgccgctgctaatcgttggcacccttcgcgctgtctcgtgcttctgctaaagtctctagccatttaaaattttgataaaccagctgcgacg

1001 cttttttctggcgagatagttcttgaatgcgggccaggatctgcacactggatatttcgggttttggggccggggcgaggggcccgtgctccc

1101 agcgccacatgcttcggcgaggcggggcctgcgagcgcggccaccgagaatcggacgggggttagtctcaaactggccggcctgctctggtgctggcctcgc

1201 gccgcccgtgatcgcggccctggggcgcaaggtggcccggctcgcgtagcggaaagatggccgcttcccggcctgctgcaggggagc

SmaI (1324)

1301 tcaaatggaggacgcccggcgggagagcgggagggtgagtcacccacacaaaggaaaggcctttccttccctcatccgtcgttcatgtgactcca

1401 cggagtaccgggcccgtccaggcaccctcgattagttgtcgagcttttggagtacgtcgtcttttagttggggggagggttttatgcgatggagtttcc

DraIII (1510)

1501 ccacactgagtggtgggagactgaagagttaggccagcttggcacttgcgtgtaattctccttgggaatttgcctttttagtttggatcttgcctcattc

AgeI (1673)

1601 tcaagcctcagacagtggttcaaagttttttcttccatttcagGTGTCGTGAAAACCTACCCCTAAAAGCCACCGGTAGGAGGCCAGCATGCTACGAGC

1701 TCTTTGGCTCTTCTGGATCTTGGTGGCCATAACAGTCTCTTACAGAAACGCTGTTCTGCTCAGGAGTCTCTGTATGTATGCTTCTGGGGTGTGTGAT

4▶ aLeuTrpLeuPheTrpI leLeuValAlaI leThrValLeuPheSerLysArgCysSerAlaGlnGluSerLeuSerCysAspAlaSerGlyValCysAsp

BstEII (1898)

1801 GGCCGCTCCAGGTCTTTCACCTCTATTCCCTCCGGACTCACAGCAGCCATGAAAAGCCTTGACCTGTCTTTCAACAAGATCACCTACATTGGCCATGGTG

38▶ GlyArgSerArgSerPheThrSerI leProSerGlyLeuThrAlaAlaMetLysSerLeuAspLeuSerPheAsnLysI leThrTyrI leGlyHisGlyA

1901 ACCTCCGAGCGTGTGCGAACCTCCAGTTCTGATTTTGAAGTCCAGCAGAATCAATACAATAGAGGGAGACGCTTTTATTCTCTGGCAGTCTTGAACA

71▶ spLeuArgAlaCysAlaAsnLeuGlnValLeuI leLeuLysSerArgI leAsnThrI leGluGlyAspAlaPheTyrSerLeuGlySerLeuGluHi

2001 TTTGGATTGTCTGATAATCACCTATCTAGTTTATCTCTCTCTGGGCCCTTTCTCTTTGAAATACTTAAACTAAATGGGAAATCCTTACCAG

104▶ sLeuAspLeuSerAspAsnHisLeuSerSerLeuSerSerTrpPheGlyProLeuSerSerLeuLysTyrLeuAsnLeuMetGlyAsnProTyrGln

Bsu36I (2150)

2101 ACATGGGGTAAACATCGCTTTTCCCAATCTCACAATTTACAACCCCTCAGGATAGGAAATGTAGAGACTTTCAGTGAGATAAGGAGAATAGATTTTG

138▶ ThrLeuGlyValThrSerLeuPheProAsnLeuThrAsnLeuGlnThrLeuArgI leGlyAsnValGluThrPheSerGluI leArgArgI leAspPheA

2201 CTGGGCTGACTTCTCTCAATGAACCTTGAATTAAGGCATTAAGTCTCCGAATTATCAGTCCCAAAGTCTAAAGTCATCCGGACATCCATCACCTGAC

171▶ IaGlyLeuThrSerLeuAsnGluLeuGluI leLysAlaLeuSerLeuArgAsnTyrGlnSerGlnSerLeuLysSerI leArgAspI leHisHisLeuTh

2301 TCTTCACTAAGCGAGTCTGCTTTCTGCTGGAGATTTTGCAGATATTCTGAGTCTGTGAGATATTTAGAACTAAGAGATACTAAGTGGCCAGGTTTC

204▶ rLeuHisLeuSerGluSerAlaPheLeuLeuGluI lePheAlaAspI leLeuSerSerValArgTyrLeuGluLeuArgAspThrAsnLeuAlaArgPhe

2401 CAGTTTTCACCACTGCCGTAGATGAAGTCAGCTCACCGATGAAGAAGTGGCATTCGGAGGCTCGGTTCTCACTGATGAAAGCTTTAACGAGCTCTGA

238▶ GlnPheSerProLeuProValAspGluValSerSerProMetLysLysLeuAlaPheArgGlySerValLeuThrAspGluSerPheAsnGluLeuLeuL

2501 AGCTGTTGCGTTACATCTTGAAGTGTGCGAGGTAGAGTTCGACGACTGTACCCTCAATGGGCTCGCGGATTTCAACCCCTCGGAGTCAGACGTAGTGAG

271▶ ysLeuLeuArgTyrI leLeuGluLeuSerGluValGluPheAspAspCysThrLeuAsnGlyLeuGlyAspPheAsnProSerLeuSerAspValValSe

2601 CGAGCTGGGTAAGTAGAAACAGTCACTATCCGGAGGTTGCATATCCCCAGTTCATTTGTTTTATGACCTGAGTACTGTCTATTCCTCTCGGAGAAG

304▶ rGluLeuGlyLysValGluThrValThrI leArgArgLeuHisI leProGlnPheTyrLeuPheTyrAspLeuSerThrValTyrSerLeuLeuGluLys

2701 GTGAAGCGAATCACAGTAGAGAACAGCAAGTCTTCTGGTTCCTGCTGCTTCCCAGCATTTAAAATCATTAGAATTTTACAGCTCAGCGAAAAATC

338▶ ValLysArgI leThrValGluAsnSerLysValPheLeuValProCysThrLeuAsnGlyLeuGlyAspPheAsnProSerLeuGluSerLeuAsnL

2801 TGATGGTTGAAGAATATTTGAAGAACTCAGCCTGTAAGGGAGCTGGCTTCTCTACAAACCTTAGTTTTGAGCCAGAATCATTTGAGATCAATGCAAAA

371▶ euMetValGluGluTyrLeuLysAsnSerAlaCysLysGlyAlaTrpProSerLeuGlnThrLeuValLeuSerGlnAsnHisLeuArgSerMetGlnLys

2901 AACAGGAGATTTTGTGACTCTGAAAAACCTGACCTCCCTTGACATCAGCAGGAACACTTTTCATCCGATGCCGACAGCTGTCAGTGGCCAGAAAAG

404▶ sThrGlyGluI leLeuLeuThrLeuLysAsnLeuThrLeuAspI leSerArgAsnThrPheHisProMetProAspSerCysGlnTrpProGluLys

BamHI (3030)

3001 ATGCGCTTCTGAAATTTGTCAGTACAGGGATCCGGGTGGTAAAAACGTCATTCTCAGACGCTGGAGGTGTTGGATGTTAGTAAACAACATCTTGACT

438▶ MetArgPheLeuAsnLeuSerSerThrGlyI leArgValValLysThrCysI leProGlnThrLeuGluValLeuAspValSerAsnAsnLeuAspS

BspHI (1898)

3101 CATTTCCTTGTCTTCTGCTCGGCTGCAAGAGCTCTATATTTCCAGAATAAGCTGAAAACACTCCCAGATGCTTCGTTGTTCCCTGTGTTGCTGGTCAT

471▶ erPheSerLeuPheLeuProArgLeuGlnGluLeuTyrI leSerArgAsnLeuLysThrLeuProAspAlaSerLeuPheProValLeuValMe

3201 GAAAATCAGAGAGAATGCAGTAAGTACTTCTCTAAAGACCACTGGTTCTTTCCCAAACCTGGAGACTTGGAAAGCAGCGACACCACCTTTGTTGTC

504▶ tLysI leArgGluAsnAlaValSerThrPheSerLysAspGlnLeuGlySerPheProLysLeuGluThrLeuGluAlaGlyAspAsnHisPheValCys

3301 TCCTGCGAATCTATCCTTTACTATGGAGACGCCAGCTCTGGCTCAAATCTGGTTGACTGGCCAGACAGTACCTGTGTGACTCTCCGCTCCGCTGC

538▶ SerCysGluLeuLeuSerPheThrMetGluThrProAlaLeuAlaGlnI leLeuValAspTrpProAspSerTyrLeuCysAspSerProProArgLeuH

3401 ACGGCCACAGGCTTCAGATGCGCGCCCTCGTCTGGATGTCACAGGCTGACTGGTGTCTGGAGTCTGCTGCGCTTCTCCTGTTGATCTTGCT

571▶ isGlyHisArgLeuGlnAspAlaArgProSerValLeuAlaCysHisGlnAlaAlaLeuValSerGlyValCysCysAlaLeuLeuSerLeuLeuLeu

3501 CGTAGGTGCCCTGTGCCACCATTTCACCGACTGTGGTACCTGAGAATGATGTGGGCGTGGCTCCAGGCCAAGAGGAGCCCAAGAAAGCTCCCTGCAGG  
604▶ uValGlyAlaLeuCysHisHisPheHisGlyLeuTrpTyrLeuArgMetMetTrpAlaTrpLeuGlnAlaLysArgLysProLysLysAlaProCysArg  
3601 GACGTTTGCATGATGCCTTTGTTTCCCTACAGTGAGCAGGATTCCCATTTGGGTGGAGAACCCTCATGGTCCAGCAGCTGGAGAACCCTGACCCGCCCTTTA  
638▶ AspValCysTyrAspAlaPheValSerTyrSerGluGlnAspSerHisTrpValGluAsnLeuMetValGlnGlnLeuGluAsnSerAspProProPheL  
ClaI (3759)  
3701 AGCTGTGTCTCCACAAGCGGGACTTCGTTCCGGGCAAAATGGATCATTGACAACATCATCGATTCCATCGAAAAAGGCCACAAAACCTGTGTTCTGCTTTTC  
671▶ ysLeuCysLeuHisLysArgAspPheValProGlyLysTrpIleIleAspAsnIleIleAspSerIleGluLysSerHisLysThrValPheValLeuSe  
3801 TGAGAACTTCGTACGGAGCGAGTGGTCAAGTACGAACTGGACTTCTCCACTTCAGGCTCTTTGACGAGAACAACGACGCGGCCATCCTTGTGTTTCTGCTG  
704▶ rGluAsnPheValArgSerGluTrpCysLysTyrGluLeuAspPheSerHisPheArgLeuPheAspGluAsnAsnAspAlaAlaIleLeuValLeuLeu  
Eco47III (3932) XcmI (3970)  
3901 GAGCCCATGAGAGGAAAGCCATTCCCCAGCGCTTCTGCAAACTGCGCAAGATAATGAACACCAAGACCTACCTGGAGTGGCCCTTGGATGAAGGCCAGC  
738▶ GluProIleGluArgLysAlaIleProGlnArgPheCysLysLeuArgLysIleMetAsnThrLysThrTyrLeuGluTrpProLeuAspGluGlyGlnG  
AvrII (4041) NheI (4066)  
4001 AGGAAGTGTGTTTGGGTAATCTGAGAAGTCAATAAAGTCTAGGTTCTCCACCCAGTTCCTGAGCTAGCTGGCCAGACATGATAAGATACATTGATGAG  
771▶ InGluValPheTrpValAsnLeuArgThrAlaIleLysSer...  
4101 TTTGGACAAACCACAAC TAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAAC  
MfeI (4215)  
4201 AAGTTAACAACAACATTCATTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGA  
PacI (4310)  
4301 AATGTTAATTAAGTCCATGACCAAAATCCCTTAACGTGAGTTTTTCGTCCACTGAGCGTCAGACCCCGTAGAAAAGATCAAAGGATCTTCTTGAGATC  
4401 CTTTTTTTTCTGCGGTAATCTGCTGTTGCAACAAAAAACACCGCTACCAGCGTGGTTTGTTCGCGGATCAAGAGTACCAACTCTTTTTCCGAA  
4501 GGTAAGTGGTTCAGCAGAGCGCAGATACCAATACTGTTCTTCTAGTGTAGCCGTAGTTAGGCCACCCTCAAGAACTCTGTAGCACCCTACATAC  
4601 CTCGCTCTGTAATCTCTGTACCAGTGGTCTGCCAGTGGGATAAGTCTGTCTTACCGGTTGGACTCAAGACGATAGTTACCGGATAAGGCGCAGC  
ApaI (4724)  
4701 GGTCCGGCTGAACGGGGGTTCTGTCACACAGCCAGCTTGGAGCGAACGACCTACACCGAACTGAGATACCTACAGCGTGAAGTATGAGAAAGCGCCAC  
4801 GCTTCCCGAAGGGAGAAAGCGGCAGGATCCCGTAAGCGGCAGGGTCCGAAACAGGAGCGCACGAGGGAGCTTCCAGGGGAAACCGCTGGTATCTT  
4901 TATAGTCTGTGCGGTTTCGCCACCTCTGACTTGAAGCGTGGATTTTTGTGATGCTCGTCAGGGGGCGGAGCCTATGGAAAAACGCCAGCAACCGCGCT  
PacI (5050)  
5001 TTTTACGGTTCCTGGCCTTTTGTGCTGCTTTTGTCTCACATGTTCTTAATTAACCTGCAGGGCCTGAAATAACCTCTGAAAGAGAACTTGGTTAGTACC  
5101 TTCTGAGGCTGAAAGAACCAGCTGTGGAATGTGTGTCACTAGGTTGGTGGAAAGTCCCAGGCTCCCAGCAGGCAGAAAGTATGCAAAGCATGCATCTCA  
5201 ATTAGTCAGCAACCAGTGTGGAAAGTCCCAGGCTCCCAGCAGGCAGAAAGTATGCAAAGCATGCATCTCAATTAGTCAGCAACCATAGTCCCAGTACT  
NotI (5329) SmaI (5358)  
5301 TCCGCCAGAGCGCGGAGGGCCTCCAGCGCGGCCCTCCCCACAGCAGGGCGGGTCCCAGCGCCACCAGGAGGCGGGCTCGGGCGGGCGGGCGC  
5401 TGATTGGCCGGGGCGGCTGACGCCGACGGCTATAAGAGACCACAAGCGACCCGAGGGCCAGACGTTCTTCGCCGAAGCTTCCGCTCAGAACGCAG  
5501 gtgaggggcggggtgtggcttccgcgggcccgcgagctggaggtcctgctccgagcggggccgggccccgctgtcgtcggcggggattagctgcgagcattc  
5601 ccgcttcgagttgcgggcgggcggggagggagagtgcgagggcctagcggcaacccctagcctcgctcgtgctccgcttgaggcctagcgtggtgtccg  
5701 cgcccgcccgctgctactccggcgcactctggtctttttttttttgttgtgttgcctgctgcctcgattgccgttcagcaataggggctaaaca  
5801 aaggagggtgcggggcttgcctcgcccggagccggagaggtcatggttggggaggaatggagggacaggagtgccggctggggccccgccccttcgga  
5901 gcacatgtccgacgccacctggatggggcgaggctggggtttttcccgaagcaaccagctggggttagcgtgccgagggcattgtggccccagcaccg  
SmaI (6102)  
6001 gcacgatctggcttggcgggcggcgttggcctgcctccctaaactagggtgaggccatcccgtccggcaccagttgcgtgctggaaagatggccgctcc  
6101 cgggcccgtgttcaaggagctcaaatggaggacggcgagcccgggtggagcgggcccgggtgagtcacccacaaaaggaagaggcctggtccctcaccg  
6201 gctgctgcttccctgtgaccccgtggtcctatcgggccgaatagtcacctcgggcttttgagcagggctagtcgcccggggggaggggatgtaatggcgt  
XcmI (6348)  
6301 tggagtttgttcacatttgggtgggtggagactagtcaggccagcctggcgtggaagtcattttggaaatttgcctccttgagtttgagcggagctaat  
BspHI (6482)  
6401 tctcgggcttcttagcggttcaaaggtatcttttaaaccttttttagGTGTTGTGAAAACCACCCTAATTCAAAGCAATCATGACTAAACCAAAATTC  
▶ MetThrLysProAsnSer  
6501 CTCATCTTCTACTGTATCATTGTTTTAGGACTGACACTTATGAAAATCCAATTATCTGAGAAATGTGAGCTTATCATAAAGAGGCCAAACGCAACCTTA  
7▶ LeuIlePheTyrCysIleIleValLeuGlyLeuThrLeuMetLysIleGlnLeuSerGluLysCysGluLeuIleIleLysArgProAsnAlaAsnLeuT  
BglII (6642)  
6601 CCAGAGTGCACAAGGACCTACCTTGAACAACACTACTTTAGATCTATCACAACAATATATCTGAGCTTACAGACTTCTGACATCCTCTCATTGTCCAA  
40▶ hrArgValProLysAspLeuProLeuGlnThrThrLeuAspLeuSerGlnAsnAsnIleSerGluLeuGlnThrSerAspIleLeuSerLeuSerLys  
6701 GCTGAGGTCCTGATAATGTCCTACAACAGACTCCAGTATCTTAATATCAGTGTTCAAATTCACACAGAGCTGGAATATTTGGATTGTCCCAAT  
73▶ sLeuArgValLeuIleMetSerTyrAsnArgLeuGlnTyrLeuAsnIleSerValPheLysPheAsnThrGluLeuGluTyrLeuAspLeuSerHisAsn  
6801 GAGCTAAAGGTGATCTTGTGCCACCAACAGTCAAGCCTCAAGCATTGGACCTCCTTTAATGCCTTTGATGCCCTGCTATATGCAAGAATTTGGCA  
107▶ GluLeuLysValIleLeuLysHisProThrValSerLeuLysHisLeuAspLeuSerPheAsnAlaPheAspAlaLeuProIleCysLysGluPheGlyA  
6901 ACATGTCCTCAACTACAGTTCCTGGGTTGAGCGGTTCTCGGTACAAAAGTTCAAAGTGTGACGCTGATTGCTCATTGAACATCAGTAAGGTTTTGCTGTT  
140▶ snMetSerGlnLeuGlnPheLeuGlyLeuSerGlySerArgValGlnSerSerValGlnLeuIleAlaHisLeuAsnIleSerLysValLeuLeuVa

7001 GTTAGGAGATGCTTATGGGAAAAAGAACCCCGAATCTCTTCGGCACGTTAGACTGAGACTCTGCATATTGTTTTCCCGTCGAAAAAGAAATCCGGT  
173▶ ILeuGlyAspAlaTyrGlyGluLysGluAspProGluSerLeuArgHisValSerThrGluThrLeuHisI leValPheProSerLysArgGluPheArg  
PvuI (7133)  
7101 TTTCTTCTGGATGTGTCGCTCAGCACTACGATCGGTTTGGAACTGTCTAACATCAAGTGTGTGCTTGAAGACCAGGGCTGCTCTTATTTCTTACGTGCTT  
207▶ PheLeuLeuAspValSerValSerThrThrl leGlyLeuGluLeuSerAsnI leLysCysValLeuGluAspGlnGlyCysSerTyrPheLeuArgAlaL  
7201 TGTCAAAGCTTGGAAAGAATCTGAAGCTCTCAAATCTTACCTGAACAATGTGGAAACAACGTGGAATTCCTTCATTAATATCCTCCAGATAGTTTGGCA  
240▶ euSerLysLeuGlyLysAsnLeuLysLeuSerAsnLeuThrLeuAsnAsnValGluThrThrTrpAsnSerPheI leAsnI leLeuGlnI leValTrpHi  
7301 TACGCCAGTCAAATATTTCTCAATTTCAAATGTGAAGCTACAAGGTCAACTTGCCTCAGGATGTTCATTATTCTGACACTTCTGAGGCTTTGTCG  
273▶ sThrProValLysTyrPheSerI leSerAsnValLysLeuGlnGlyGlnLeuAlaPheArgMetPheAsnTyrSerAspThrSerLeuLysAlaLeuSer  
7401 ATACATCAAGTTGTCACGTGTCTTCAGCTTCCCCAAAGTTACATATACAGTATCTTTGCCAATATGAACATCCAAAACCTTACAATGTCTGGAACAC  
307▶ I leHisGlnValValThrAspValPheSerPheProGlnSerTyrI leTyrSerI lePheAlaAsnMetAsnI leGlnAsnPheThrMetSerGlyThrH  
7501 ACATGGTCCACATGCTGTGCCCGTCCCAAGTTAGCCATTTCTGCATGTGGACTTTACAGATAACCTTTAAACAGACATGGTTTTTAAAGACTGTAGAAA  
340▶ isMetValHisMetLeuCysProSerGlnValSerProPheLeuHisValAspPheThrAspAsnLeuLeuThrAspMetValPheLysAspCysArgAs  
7601 CTTAGTTAGATTGAAAACTTAGTTTACAAAAAGAAATCAGTTAAAAAACCTTGAGAAATATAATCCTCACATCTGCAAGATGACATCCCTACAAAACTA  
373▶ nLeuValArgLeuLysThrLeuSerLeuGlnLysAsnGlnLeuLysAsnLeuGluAsnI leI leLeuThrSerAlaLysMetThrSerLeuGlnLysLeu  
BstBI (7786)  
7701 GACATTAGCCAGAATCTCTAAGGTACAGCGATGGGGGAATCCCATGCGCTGGACCAGAGTTTGTAGTTTTAAATTTGTCTTCGAATATGCTTACAG  
407▶ AspI leSerGlnAsnSerLeuArgTyrSerAspGlyGlyI leProCysAlaTrpThrGlnSerLeuLeuValLeuAsnLeuSerSerAsnMetLeuThrG  
7801 GCTCTGTCTTCAGATGCTTACTCCCAAAGTCAAGTCTTGCACCTTCAACAACAAGGATAATGAGCATCCCTAAAGATGCACCCACTGCAGGCTTT  
440▶ lySerProSerArgCysLeuProProLysValLysValLeuAspLeuThrLeuGlyHisAsnAsnArgI leMetSerI leProLysThrCysValTheHisLeuAlaLe  
7901 GCAGGAACCTCAATGTAGCATCAACTCCTTAAGTACCTTCTGGTGGCGGGCCTTACAGCCTTTCTGTGCTGGTCATCGACCATAACTCAGTTTCC  
473▶ uGlnGluLeuAsnValAlaSerAsnSerLeuThrAspLeuProGlyCysGlyAlaPheSerSerLeuSerValLeuVal I leAspHisAsnSerValSer  
8001 CATCCCTCTGAGGATTTCTCCAGAGCTGTCAGAATATTAGATCCCTAACAGCGGGAACAACCCATTCCAATGCACATGTGAGCTGAGGGACTTTGTCA  
507▶ HisProSerGluAspPhePheGlnSerCysGlnAsnI leArgSerLeuThrAlaGlyAsnAsnProPheGlnCysThrCysGluLeuArgAspPheValI  
8101 AGAACATAGGCTGGGTAGCAAGAGAAGTGGTGGAGGGCTGGCTGACTTTACAGTGTGACTACCCAGAAAGCTCTAGGGGAAGTGCAGTGGGGACTT  
540▶ ysAsnI leGlyTrpValAlaArgGluValValGluGlyTrpProAspSerTyrArgCysAspTyrProGluSerSerArgGlyThrAlaLeuArgAspPh  
8201 CCACATGTCTCCACTATCCTGTGATACTGTTCTGTGACTGTCACCATCGGGCCACTATGCTGGTGTGGTGTGACTGGGGCTTCTCTGTCTCTAC  
573▶ eHisMetSerProLysCysAspThrValLeuLeuThrValLeuLeuAlaValI leMetSerI leProLysThrCysValTheHisLeuAlaLeuTrp  
8301 TTTGACCTGCCCTGGTATGTGAGGATGCTGTGTCAGTGGACACAGACCAGGCACAGGGCCAGGCACATCCCTTAGAGGAAGTCCAGAGAAACCTCCAGT  
607▶ PheAspLeuProTrpTyrValArgMetLeuCysGlnTrpThrGlnThrArgHisArgAlaArgHisI leProLeuGluGluLeuGlnArgAsnLeuGlnP  
8401 TCCATGCTTTTGTCTCATACTGAGTGGCATGATTCTGCCTGGTGAAGAACAATTACTACCAACCTAGAGAAAGATGACATCCAGATTTGCCCTCCATGA  
640▶ heHisAlaPheValSerTyrSerGlyHisAspSerAlaTrpValLysLeuThrValLeuThrI leGlyHisAsnAsnArgI leMetSerI leProLysThrCysValTheHisG  
8501 GAGGAACCTTTGTCCTGGCAAGAGCATTGTGGAGAACATCAATTTTCATTGAGAAGAGTTACAAGTCCATCTTTGTGCTGTCTCCCACTTCCATCCAG  
673▶ uArgAsnPheValProGlyLysSerI leValGluAsnI leI leAsnPheI leGluLysSerTyrLysSerI lePheValLeuSerProHisPheI leGln  
8601 AGTGAGTGGTGCATTATGAAGTCTATTTTGGCCATCAATCTCTCCATGAAGGCTCTGATAACTTAATCTCATCTTTGCTGGCACCATTCCCGAGT  
707▶ SerGluTrpCysHisTyrGluLeuTyrPheAlaHisHisAsnLeuThrValLeuThrI leGlyHisAsnAsnArgI leMetSerI leProLysThrCysValTheHisG  
8701 ACTCCATCCCTACCAATTACCACAAGCTCAAACTCTCATGTCACGAAGGACCTATCTGGAATGGCCACAGAGAAGAACAAGCATGGACTTTTTTGGGC  
740▶ yrSerI leProThrAsnTyrHisLysLeuLysThrLeuMetSerArgArgThrTyrLeuGluTrpProThrGluLysAsnLysHisGlyLeuPheTrpAl  
8801 AAACCTAAGAGCATCCATTAATGTAAAGTGGTAAACAGGCAGAAAGGAACTGTTACACACAGCAATAAGAATATCCACCGCTAGGAGCAGGTTTCCCG  
773▶ aAsnLeuArgAlaSerI leAsnValLysLeuValAsnGlnAlaGluGlyThrGlnGln • • •  
XbaI (8948)  
8901 AATGACACAAAACGTGCAACTTGAAGTCCGCCTGGTCTTTCCAGGTCTAGAGGGGTAACACTTTGTACTGCGTTTGGCTCCACGCTCGATCCACTGGCG  
9001 AGTGTTAGTAACAGCACTGTTGCTTCTGAGCGGAGCATGACGGCCGTGGGAACCTCCTCTTGGTAAACAAGGACCCACGGGGCCAAAAGCCACGCCCACAC  
9101 GGGCCCGTCATGTTGCAACCCAGCACGGCGACTTTACTGCGAAACCACTTTAAAGTGACATTGAAACTGGTACCACACACTGGTGACAGGCTAAGG  
Eco47III (9274)  
9201 ATGCCCTTCAGTACCCGAGGTAACACGGCACTCGGGATCTGAGAAGGGGACTGGGGCTTCTATAAAAGCGCTCGGTTAAAAAGCTTCTATGCCTG  
BstEII (9306)  
9301 AATAGGTGACGGAGGTGCGCACCTTTCTTTGCAATTACTGACCCTATGAATACACTGACTGTTTGACAATTAATCATCGGCATAGTATATCGGCATAG  
9401 TATAATACGACTCACTATAGAGGGCCACCATGAAGACCTTCAACATCTCTCAGCAGGATCTGGAGCTGGTGGAGGTGCCACTGAGAAGATCACCATGC  
1▶ MetLysThrPheAsnI leSerGlnGlnAspLeuGluLeuValGluValAlaThrGluLysI leThrMetL  
9501 TCTATGAGGACAACAAGCACCATTGTCGGGGCGGCCATCAGGACCAAGACTGGGGAGATCATCTCTGCTGTCCACATTGAGGCTACATTGGCAGGGTCCAC  
24▶ euTyrGluAspAsnLysHisHisValGlyAlaAlaI leArgThrLysThrGlyGluI leI leSerAlaValHisI leGluAlaTyrI leGlyArgValTh  
9601 TGTCTGTGCTGAAGCCATTGCCATTGGGCTGTGCTGTGAGCAACGGGCGAAGGACTTTGACACCATTGTGGCTGTGAGGCACCCCTACTCTGATGAGGTG  
57▶ rValCysAlaGluAlaI leAlaI leGlySerAlaValSerAsnGlyGlnLysAspPheAspThrI leValAlaValArgHisProTyrSerAspGluVal  
9701 GACAGATCCATCAGGGTGGTCCAGCCCTGTGGCATGTGCAGAGAGCTCATCTGACTATGCTCCTGACTGCTTTGTGCTCATTGAGATGAATGGCAAGC  
91▶ AspArgSerI leArgValValSerProCysGlyMetCysArgGluLeuI leSerAspTyrAlaProAspCysPheValLeuI leGluMetAsnGlyLysL  
9801 TGGTCAAAACCACTTGGAGAACTCATCCCCCTCAAGTACACCAGGAACCTGAATTAATTCGCTAGGATTATCCCTAATACCTGCCACCCCACT  
124▶ euValLysThrThrl leGluGluLeuI leProLeuLysTyrThrArgAsn • • •  
MfeI (9943)  
9901 CTTAATCAGTGGTGAAGAACGGTCTCAGAAGTGGTTTTCATTTGGCCATTAAGTTTAGTAGTAAAAGACTGGTTAATGATAACAATGCATCGTAAA  
10001 ACCTTCAGAAGAAAGGAGAATGTTTTGTGACCACCTTGGTTTTCTTTTTTGGCTGTGGCAGTTTAAAGTTATTAGTTTTTAAATCAGTACTTTTTAA  
10101 TGGAACAACCTTGACAAAAATTTGTACAGAATTTGAGACCCATTAAGAAAGTAAATGAGAAACCTGTGTCTCTTTGGTCAACCCGAGACATTT  
10201 AGGTGAAAGACATCTAATCTGGTTTTACGAATCTGGAAACTTCTTGAAGTAAATCTTGAGTTAACACTTCTGGGTGAGAAATAGGGTTGTTTTCC  
10301 CCCACATAATTGGAAGGGGAAGGAATATCATTTAAAGCTATGGGAGGTTTCTTTGATTACAACACTGGAGAGAAATGACAGCATGTTGCTGATTGCCTGT  
10401 CACTAAAACAGGCCAAAACTGAGTCTTGGTTCATAGAAGCTG