

pTRIOZ-hIgG1

Plasmid for high yield production of recombinant human IgG1 kappa mAbs

Catalog code: ptrioz-higg1

<https://www.invivogen.com/ptrioz-higg1>

For research use only

Version 21E28-ED

PRODUCT INFORMATION

Contents

- 20 µg of pTRIOZ-hIgG1 plasmid provided as lyophilized DNA
- 1 ml of Zeocin™ (100 mg/ml)

Storage and Stability

- pTRIOZ-hIgG1 is provided as a lyophilized powder and shipped at room temperature. Upon receipt, store product at -20°C.
- Store resuspended product at -20°C. Resuspended product is stable for at least 1 year when properly stored.
- Avoid repeated freeze-thaw cycles.
- Store Zeocin™ 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.

PRODUCT DESCRIPTION

The pTRIOZ plasmid collection has been designed specifically for high yield production of whole recombinant monoclonal antibodies (mAbs).

The pTRIOZ plasmids contain three distinct cassettes for the expression of the heavy and light chain of the mAb as well as antibiotic selection with Zeocin™ in both bacterial (such as *E. coli*) and mammalian (such as CHO) cells. Each cassette is under the control of unique composite promoters for optimal expression (see *Plasmid features for more details*). For successful mAb production, a precise expression ratio of the heavy to light chain is required¹. In the pTRIOZ plasmids this important ratio is under the control of the human ferritin heavy (FerH) and light (FerL) chain promoters, which natively drive the successful co-expression of the two ferritin subunits². Additionally, the pTRIOZ plasmids contain unique multiple cloning sites (MCS) upstream of both the heavy and light chain constant (CH and CL) regions. This enables the cloning of variable (VH and VL) regions of any given antibody.

Majority of mAbs are produced by recombinant DNA technology in mammalian cells, either through transient or stable gene expression. The pTRIOZ plasmid collection has been designed to be used for either method. Transient or stable transfection of mammalian cell lines, such as CHO cells, with a recombinant pTRIOZ plasmid results in high-yield production of an IgG mAb that can be purified from the supernatant using an appropriate Protein A or Protein G affinity chromatography method.

pTRIOZ-hIgG1 expresses the constant region of the heavy (CH) chain from human IgG1 and the constant region of the human kappa light chain (CL). pTRIOZ-hIgG1 is selectable in both bacterial and mammalian cells with Zeocin™.

PLASMID FEATURES

CASSETTE 1: mAb HEAVY CHAIN

- **AldA enh/ hFerH:** This composite promoter combines the human aldehyde dehydrogenase (aldA) enhancer and the core promoter of the human ferritin heavy chain gene.
- **MCS1:** To facilitate cloning of the variable heavy (VH) chain, the multiple cloning site contains the following restriction sites that are compatible with many different enzymes, 5'- *AgeI, MluI, EcoRV, and NheI* -3'.
- **hIgG1:** The constant region of the human immunoglobulin IgG1 heavy chain.
- **βGlo pAn:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription.

CASSETTE 2: mAb LIGHT CHAIN

- **hCMV enh / hFerL prom:** This composite promoter combines the human cytomegalovirus (CMV) immediate-early gene 1 enhancer and the core promoter of the human ferritin light chain gene.
- **MCS2:** To facilitate cloning of the variable light (VL) chain, the multiple cloning site contains the following restriction sites that are compatible with many different enzymes, 5'- *SgrAI, AscI, PmeI, NcoI, and BsiWI* -3'.
- **Human κ light chain:** The constant region of the human kappa light chain
- **SV40 pAn:** The Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA

CASSETTE 3: Zeocin™ SELECTION

- **mCMV/hEF1-HTLV prom:** This composite promoter combines mouse cytomegalovirus (mCMV) immediate-early gene 1 enhancer, the elongation Factor-1α (EF-1α) core promoter, as well as the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) type 1 long terminal repeat. The EF-1α promoter exhibits a strong activity and yields long lasting expression of a transgene *in vivo*. The R-U5' has been coupled to the EF-1α core promoter to enhance stability of RNA.
- **EM7 prom:** This is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*. EM7 is located within an intron and is spliced out in mammalian cells.
- **Sh Ble gene:** Resistance to Zeocin™ is conferred by the *Sh ble* gene from *Streptoalloteichus hindustanus*. The same gene confers resistance in both mammalian cells and *E. coli*.
- **hEF-1α pAn:** This provides a strong polyadenylation signal. InvivoGen uses a sequence that starts after the stop codon of the EF1 cDNA and finishes after a bent structure rich in GT.

GENERAL FEATURES: pTRIOZ-hIgG1

- **5' UTR:** The 5' UTR enhances mRNA stability and protein translation.
- **Ori:** A minimal *E. coli* origin of replication.

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



PLASMID RESUSPENSION

- Centrifuge the tube containing the lyophilized pTRIOZ-hlgG1 plasmid to pellet the DNA.
- To obtain a plasmid solution at 1 µg/µl, resuspend the DNA in 20 µl of sterile endotoxin-free H₂O.
- Store resuspended plasmid at -20°C.

GENERAL METHODS

Obtaining the VH and VL sequences

To obtain the cDNA sequence of the variable heavy (VH) and light (VL) regions from an antibody producing hybridoma, total RNA or mRNA is extracted and reverse-transcribed to cDNA. PCR is performed with 5' degenerate primers to anneal to the unknown VH and VL regions and the 3' primers designed to anneal to the "known" CH and CL regions. The resulting amplicons must be sequenced. Additionally, the VH and VL chains of the mAb can be commercially synthesised. This allows for codon optimization, both for the expression system, as well as ensuring that restriction sites in the MCS are avoided. Furthermore, the 5' and 3' cloning ends for both the VH and VL chain regions can be added.

Cloning mAb variable regions into pTRIOZ

Plasmid amplification and cloning can be performed in *E. coli* GT116 or other commonly used laboratory strains such as DH5α. For selection in *E. coli*, Zeocin™ is commonly used at 25 µg/ml in liquid or solid media

- Variable Heavy (VH) chain

In pTRIOZ-hlgG1, the constant region of the human IgG1 heavy chain is preceded by a MCS containing four restriction sites: AgeI, MluI, EcoRV, and NheI. We recommend using the AgeI restriction site for insertion of the 5' end of the mAb VH chain (including the native signal sequence).

In pTRIOZ-hlgG1, NheI must be used for insertion of the 3' end of the VH chain to maintain the integrity of the constant region. Therefore, we recommend to introduce an NheI site at the 3' end of the variable region, in frame with the constant region of the human IgG1 heavy chain. This ensures that no additional amino acids are introduced into the mAb sequence.

- Variable Light (VL) chain

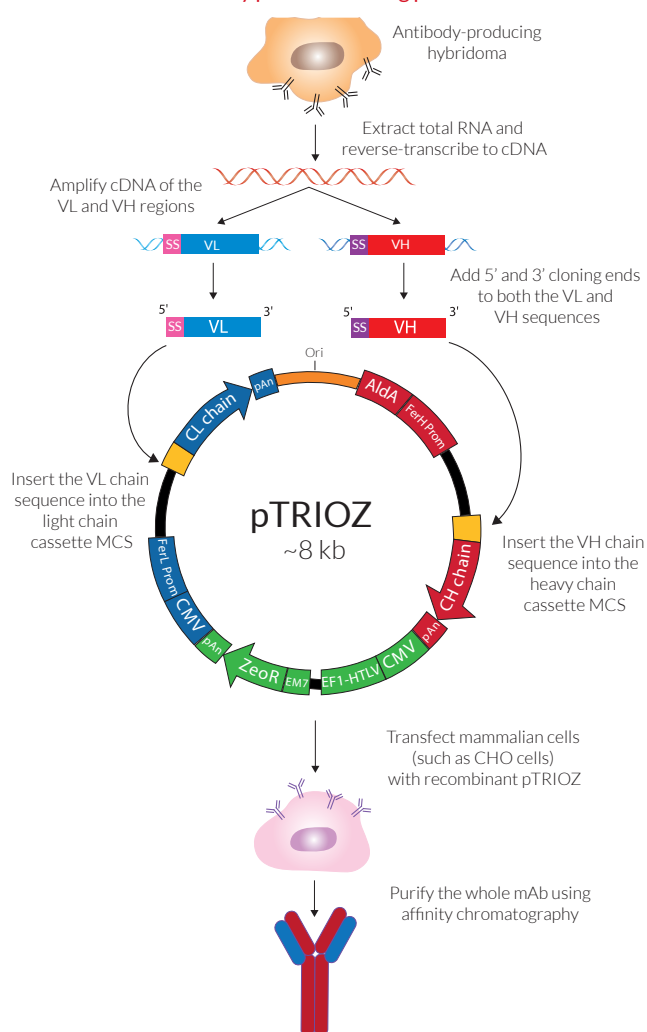
In pTRIOZ-hlgG1, the constant region of the human kappa light chain is preceded by a MCS containing five restriction sites: SgrAI, Ascl, PmeI, NcoI, and BsiWI. We recommend using the SgrAI restriction site for insertion of the 5' end of the mAb VL chain (including the native signal sequence).

In pTRIOZ-hlgG1, BsiWI must be used for insertion of the 3' end of the VL chain to maintain the integrity of the constant region. Therefore, we recommend to introduce an BsiWI site at the 3' end of the VL chain, in frame with the constant region of the human kappa light chain. This ensures that no additional amino acids are introduced into the mAb sequence.

Antibody production

The pTRIOZ plasmid collection is designed for mAb production in transient-expressing CHO and HEK cells as well as for establishing stable-expressing cell lines. Specifically for stable-expressing cell lines, 72 hours after transfection, cells should be placed into fresh medium containing 50-200 µg/ml of Zeocin™, the selection antibiotic. *Note: The optimal Zeocin™ concentration for selection should be calculated by seeding native CHO cells with different concentrations of Zeocin™ and monitoring both cell growth and viability.*

Antibody production using pTRIOZ



The selection medium should be changed every 2-3 days until cell viability and growth both become stable. Zeocin™-resistant stable cell pools are obtained typically between 7-10 days after selection. The selected stable cell pools can be used for bioproduction of mAbs in batch, fed batch or perfusion process modes.

Antibody purification

The resulting mAb can be purified from the supernatant using the appropriate Protein A or Protein G affinity chromatography.

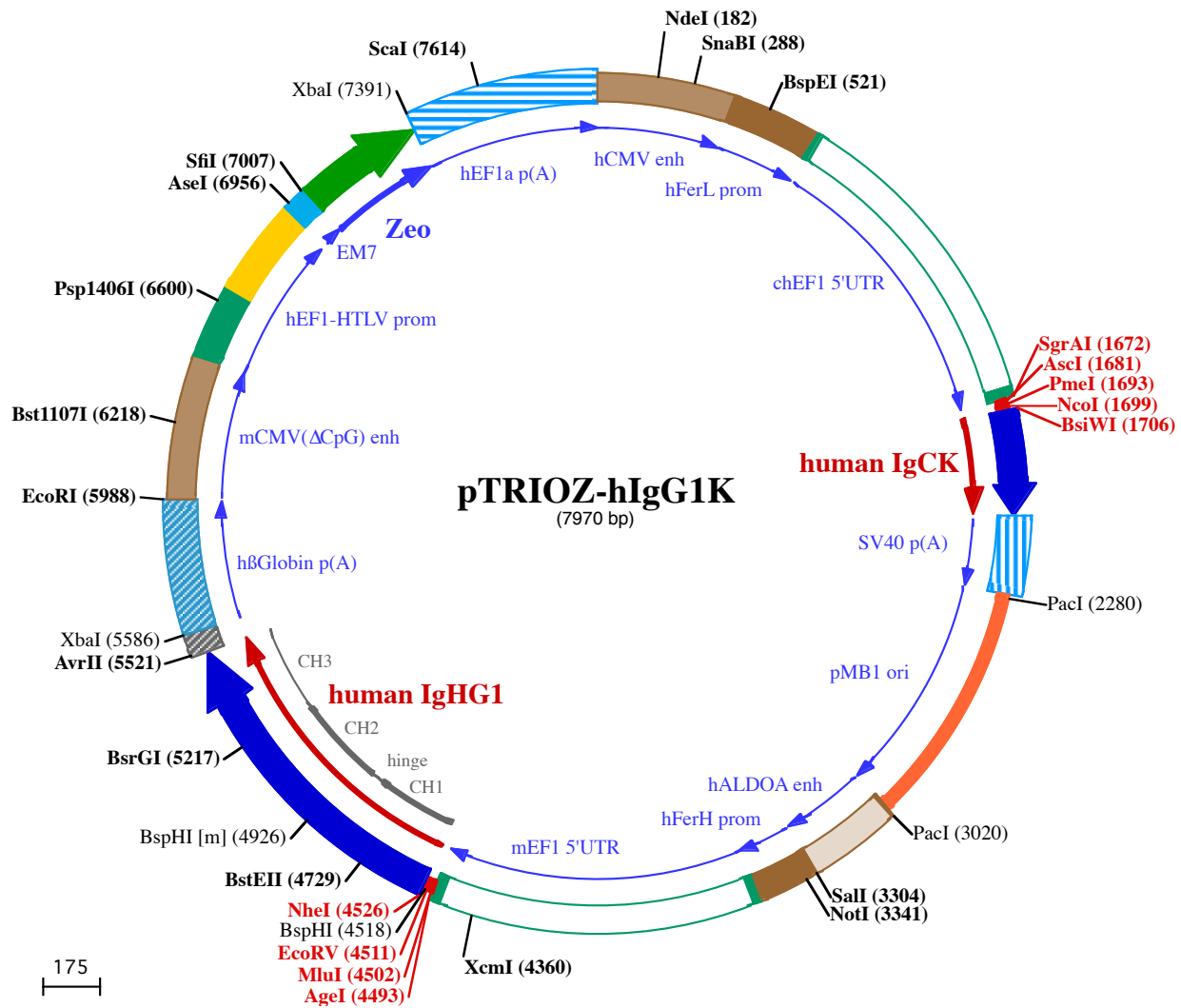
1. Prentice, H.L. et al., 2007. High level expression of proteins using sequences from the ferritin heavy chain gene locus. *J Biotech.* 128:50-60. 2. Rita costa, A. et al., 2010. Guidelines to cell engineering for monoclonal antibody production. *Eur J Pharm Biopharm.* 74(2):127-138.

RELATED PRODUCTS

Product	Catalog Code
ChemiComp GT116	gt116-11
LyoVec™	lyec-12
Protein G / Agarose	gel-agg-5
Zeocin™	ant-zn-1

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 CCTGCAGGCGTTACATAACTTACGGTAAATGGCCCGCTGGCTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAA

101 CGCCAATAGGGACTTTCCATTGACGTCAATGGGTGGAGTATTTACGGTAAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCC

201 TATTGACGTCAATGACGGTAAATGGCCCGCTGGCATTATGCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATC

301 GCTATTACCATGATGATGCGGTTTTGGCAGTACATCAATGGGCGTGGATAGCGGTTTGACTCACGGGGATTCCAAGTCTCCACCCCAATTGACGTCAATG

401 GGAGTTTGTTTTACTAGTCAGGGCCCCAACCCCCCAAGCCCCATTTACAACACGCTGGCGCTACAGGCGCGTGACTTCCCCTTGCTTTGGGGCGGG

501 GGGCTGAGACTCTATGTGCTCCGATTGGTCAGGCACGGCTTCGGCCCGCTCTGCCACCGCAGATTGGCCGCTAGGCCTCCCCGAGCGCCTGCC

601 TCCGAGGGCCGGCGCACCATAAAAGAAGCCGCCCTAGCCACGTCCCCTCGAGTTCGGCGGTCCCGCGGGTGTCTCAAGCTTGCCGCGAGAACAGG

701 taagtgccgtgtgtggttcccgcgggcctggcctctttacgggttatggccttgcgtgcctgaattacttccatgccccggctgcagtacgtgattc

801 ttgatcccagacttccgggttgaagtgggtgggagagtccagggccttgcgcttaaggagcccttcgctcgtgcttgagttgaggcctggcttggcg

901 ctggggccgcccgtgctaactctggtggcaccttcgcccctgctcgtgcttgcgtaagtctctagcatttaaaatgttataaccagctgcgagc

1001 cttttttctggcgagatagtctttaaataatgcccagatctgcacactggtatctcggtttttggggccgcccggcgagggcccgctgcgtccc

1101 agcgcacatgctcggcgaggcgggctgcgagcggccaccgagaatcggacggggtagtctcaactggccggcctgctcgtgctgctggcctgc

1201 gccgcctgtatcgcgccctgggcggcaaggctggcccggcggcaccagttcgtgagcggaaagatggccgcttcccggcctgctgcagggagc

1301 tcaaatggaggagcggcgcccgggagagcgggaggtagtcaccacacaaggaaaggccttctctcatcctcgtcgttcatgtgactcca

1401 cggagtaccggcgccgctccaggcacctcgattagttgctcagcctttggagtacgtcgtccttaggtggggggaggggtttatgcatggagttcc

1501 ccacactgagtgggtggagactgaagagttaggccagcttggcacttgatgtaattctccttgaatttgcctttttgagtttgatcttgcctcttc

1601 tcaagcctcagacagtggttcaagttttttcttccatttcagGTGTCGTGAAAACCTACCCCTAAAAGCCACCGGCGAGGCGGCCAAGTTTAAACACC

1701 ATGGAAACGTACGGTGGCTGCACCATCTGCTTTCATCTTCCCGCATCTGATGAGCAGTTGAACTCTGGAACCTGCCTCTGTTGTGTCCTGCTGAATAACT

1801 TCTATCCCAGAGAGGCCAAAGTACAGTGAAGTGGATAACGCCCTCAATCGGGTAACTCCAGGAGAGTGTACAGAGCAGGACAGCAAGGACAGCAC

1901 CTACAGCCTCAGCAGCACCCTGACGCTGAGCAAAGCAGACTACGAGAAACCAAAAGTCTACGCTGCGAAGTACCCATCAGGGCCTGAGCTCGCCGTC

2001 ACAAGAGCTTCAACAGGGGAGAGTGTAGAGGGACCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTG

2101 AAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAACATTCATTCTTT

2201 ATGTTTCAGGTTACGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATGTTAATTAAGTACCATGACCAAAT

2301 CCTTAACGTGAGTTTTCTGTTCACTGAGCGTCAGACCCGTAGAAAAGATCAAAGGATCTTCTTGTGATCCTTTTTTCTGCGGTAATCTGCTGCTTG

2401 CAAACAAAAAACACCCTACCAGCGGTGGTTTGTGGCCGATCAAGAGCTACCAACTCTTTTTCCGAAGTAACTGGCTTACGAGAGCGCAGATAC

2501 CAAATACTGTTCTTAGTGTAGCCGTAGTTAGCCACCCTCAAGAACTCTGTAGCACCGCTACATACCTCGCTCTGTAATCCTGTTACCAAGTGC

2601 TGCTGCCAGTGGCGATAAGTCGTGCTTACCGGTTGGACTCAAGACGATAGTTACCGGATAAGGCGCAGCGGTCGGGCTGAACGGGGGTTCTGTCACA

2701 CAGCCCAGCTTGAGCGAACGACCTACCCGAACCTGAGATACCTACAGCGTAGCTATGAGAAAGCGCCACGCTTCCGAAGGGAGAAAGCGGACAGGT

2801 ATCCGGTAAGCGGAGGTCGGAACAGGAGAGCGCAGGGAGCTTCCAGGGGAAACGCTGGTATCTTTATAGTCTGTGCGGTTTCGCCACCTCTG

2901 ACTTGAGCGTCGATTTTTGTGATGCTCGTCAGGGGGCGGAGCCTATGAAAAACGCCAGCAACGCGCCTTTTTACGGTTCCTGGCCTTTTGTGGCCT

3001 **Pacl (3020)**
TTTGCTCACATGTTCTTAATTAACCTGCaggcgaactcagcttccttcgtttccgacttttccatccgcgtctcacttccccgttccgcctcccc
3101 attgccaacattctggctgagtcacggcgccccagagcgccaggctgggggaaaggagcagaagggagggccctagcgaccccggggatgtggtccga
3201 gtcacgtccgaggggggtggggagggatcgtgttctcggcgcccccttcttagcgcggcctctgggctgcgcctctcggggcgccccgtagccag

3301 **SalI (3304)** **NotI (3341)**
tccgtcgACTAGTTCCGCCAGAGCGCGGAGGGCTCCAGCGGCCGCCCTCCCCACAGCAGGGGCGGGTCCC CGCICACCGGAAGGAGCGGGCTCG
3401 GGGCGGGCGCGCTGATTGGCCGGGGCGGGCTGACGCCGACGCGGTATAAGAGACCACAAGCGACCCGAGGGCCAGACGTTCTTCGCGGAAGCTTGC
3501 CGTCAGAACGCAGgtgaggggagggtgtggcttccgaggcccgagctggaggctctgctccgagcgggcccggccccgctgctgctcgggggatta
3601 gctgcgagcattccccgttcgagttgccccggcgggaggcagagtgcgaggcctagcggcaaccctgtagcctcgctcgtgctccggctgaggcct
3701 agcgtggtgtccgcgcccccgctgctactccggcgcactctggtctttttttttttttgttgttgttgcctgctgccttcgattgccgttcagca
3801 ataggggctaacaaggagggtgccccggcttgcctcgccggagcccggagaggtcatggttgggaggaatggagggacaggagtgccggctggggccc
3901 gcccgccttcggagcacatgtccgacgcacactggatggggcgaggcctggggttttccgaagcaaccaggctggggtagcgtgcgagggccatgtg
4001 gccccagacccggcacgatctggcttggcgcccgcttgcctgcctccctaactagggtgaggccatcccgccggcaccagttgcgtgcgtggaa
4101 agatggccgctccgggcccctgttgcaaggagctcaaatggaggacgcccagcccggaggcgggtagtcacccacacaaggaagagggcc
4201 tggtcctcaccggctgctgcttctgtgaccccggtctctatcggcgcaatagtcactcgggcttttgagcacggttagtcgcggcggggggaggg

4301 **XcmI (4360)**
gatgtaatggcgttggagttgtcacatttgggtgggtggagactagt caggccagcctggcgtggaagtcatttttggaatttgtccccttgagtttt

4401 **AgeI (4493)**
gagcggagctaattctcgggcttcttagcgggtcaaaggtatcttttaaaccttttttaggTGTTGTGAAAACACCGCTAATTCAAAGCAA **CCGGTCCG**

4501 **EcoRV (4511)** **NheI (4526)**
MluI (4502) **BspHI (4518)**
CACGCGTAGATATCACGTCATGAAAGCTAGCACCAAGGCCCATCGGTCTTCCCCCTGGCACCTCTCCAAGAGCACCTCTGGGGGCACAGCGGCCCTG
1▶ A S T K G P S V F P L A P S S K S T S G G T A A L

4601 GGCTGCCTGGTCAAGGACTACTTCCCCGAGCCGGTGACGGTGTCTGGAAGTCAAGCCCTGACCAGCGGCTGCACACCTTCCCGCTGTCTACAGT
26▶ G C L V K D Y F P E P V T V S W N S G A L T S G V H T F P A V L Q

4701 **BstEII (4729)**
CCTCAGGACTCTACTCCCTCAGCAGCGTGGTGACCGTGCCCTCCAGCAGCTTGGGCACCCAGACCTACATCTGCAACGTGAATCACAAGCCCAGCAACAC
59▶ S S G L Y S L S S V V T V P S S S L G T Q T Y I C N V N H K P S N T

4801 CAAGTGGACAAGAAAGTTGAGCCAAATCTTGTGACAAAAGTACACATGCCACCGTGCCAGCACCTGAACTCTGGGGGACCGTCAGTCTTCTCCTC
92▶ K V D K K V E P K S C D K T H T C P P C P A P E L L G G P S V F L

4901 **BspHI [m] (4926)**
TCCCCCAAACCAAGGACACCCTCATGATCTCCCGGACCCTGAGGTCACATGCGTGGTGGTGGACGTGAGCCACGAAGACCCTGAGGTCAAGTTCA
126▶ F P P K P K D T L M I S R T P E V T C V V V D V S H E D P E V K F

5001 ACTGGTACGTGGACGGCTGGAGGTGCATAATGCCAAGACAAAGCCGCGGGAGGAGCAGTACAACAGCACGTACCCTGGTGGTCAAGCTCCTCACCCTCT
159▶ N W Y V D G V E V H N A K T K P R E E Q Y N S T Y R V V S V L T V L

5101 GCACCAGGACTGGCTGAATGGCAAGGAGTACAAGTGAAGGTCTCCAACAAAGCCCTCCAGCCCCATCGAGAAAACCATCTCCAAGCCAAAGGGCAG
192▶ H Q D W L N G K E Y K C K V S N K A L P A P I E K T I S K A K G Q

5201 **BsrGI (5217)**
CCCCGAGAACCACAGGTGTACACCTGCCCCATCCCGGGACGAGTGACCAAGAACCAGGTCAGCCTGACCTGCCTGGTCAAAGGCTTCTATCCCAGCG
226▶ P R E P Q V Y T L P P S R D E L T K N Q V S L T C L V K G F Y P S

5301 ACATCGCGTGGAGTGGGAGGCAATGGCAGCCGAGAACTACAAGACCAGCCTCCCGTGTGGACTCCGACGGCTCCTTCTCTCTACAGCAA
259▶ D I A V E W E S N G Q P E N N Y K T T P P V L D S D G S F F L Y S K

5401 GCTCACCGTGACAAGAGCAGGTGGCAGCAGGGGAACGCTTCTCATGCTCCGTGATGCATGAGGCTCTGCACAACCACTACACGAGAAGAGCCTCTCC
292▶ L T V D K S R W Q Q G N V F S C S V M H E A L H N H Y T Q K S L S

5501 **AvrII (5521)** CTGTCTCCGGTAAATGAGTCTTAGGAGCAGGTTTCCCAATGACACAAAACGTGCAACTTAAAACCTCCGCTGGTCTTCCAGGCTAGAAAGCTCGCTT
326▶ L S P G K • XbaI (5586)

5601 TCTTGCTGTCCAATTTCTATTAAGGTTCTTTGTTCCCTAAGTCCAACACTAACTGAGGGATATTATGAAGGCCCTTGAGCATCTGGATTCTGCCTA

5701 ATAAAAACATTTATTTTCATTGCAATGATGTATTTAAATTATTTCTGAATATTTTACTAAAAAGGGAATGTGGGAGTCAAGTGCATTTAAAAACATAAAG

5801 AAATGAAGAGCTAGTTCAAACCTTGGGAAAATACACTATATCTTAAACTCCATGAAAGAAGGTGAGGCTGCAAACAGCTAATGCACATTGGCAACAGCCC

5901 **EcoRI (5988)** CTGATGCCTATGCCTTATTCATCCCTCAGAAAAGGATTCAAGTAGAGGCTTGATTTGGAGTTAAAGTTTTGCCATGCTGTATTTTAGAATTCTGCAGG
6001 AGTCAATGGGAAAACCCATTGGAGCCAAGTACACTGACTCAATAGGGACTTTCCATTGGGTTTTGCCAGTACATAAGGTCAATAGGGGGTGAAGTCAAC
6101 AGGAAAGTCCCATTGGAGCCAAGTACATTGAGTCAATAGGGACTTTCCAATGGGTTTTGCCAGTACATAAGGTCAATGGGAGGTAAGCCAATGGGTTTT

6201 **Bst1107I (6218)** TCCCACTACTGACATGTATACTAGTCAATTAGGGACTTTCCAATGGGTTTTGCCAGTACATAAGGTCAATAGGGGTGAATCAACAGGAAAAGTCCCATTG
6301 GAGCCAAGTACACTGAGTCAATAGGGACTTTCCATTGGGTTTTGCCAGTACAAAAGGTCAATAGGGGGTGAAGTCAACAGGAAAAGTCCCATTG
6401 ATACATAAGGTCAATAGGGGTGACTAGTCAGTGGGCAGAGCGCACATCGCCCGGAGAAGTTGGGGGGAGGGGTCGGCAATTGAACGGGTGCCTAGAGAA

6501 **Psp1406I (6600)** GGTGGCGCGGGTAAACTGGGAAAGTATGTCGTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCCGTGAA
6601 CGTTCTTTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCCACGCGCCGCCCTACCTGAGGCGCCCAT
6701 CCACGCCGTTGAGTTCGGTCTGCGCCTCCCGCCTGTGGTGCCTCCTGAAGTGCCTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACCGGGCC
6801 TTTGTCCGGCGCTCCCTTGGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCTGACCCTGCTTGCTCAACTCTACGTCTTTGTTTTGTTTTCTGT

6901 **AseI (6956)** TCTGCGCGTTACAGATCCAAGCTGTGACCGGCGCCTACAACAGTAGTTGACAATTAATCATCGGCATAGTATATCGGCATAGTATAATACGACTCACT

7001 **SfiI (7007)** ATAGGAGGGCCATCATGGCCAAGTTGACCAAGTCCGGTCCGGTGTCCACGCGCGGACGTCGCCGGAGCGGTCGAGTTCTGGACCGACCGGCTCGGGTT
7101 CTCCCGGACTTCGTGGAGGACACTTCGCTGGTGTGGTCCGGGACGACGTGACCCTGTTTCATCAGCGCGTCCAGGACCAGGTTGGTCCGGACAACACC
29▶ S R D F V E D D F A G V V R D D V T L F I S A V Q D Q V V P D N T
7201 CTGGCCTGGGTGTGGGTGCGCGGCTGGACGAGCTGTACGCCAGTGGTCCGAGGTCGTGTCCACGAACCTCCGGGACGCTCCGGGCGGCCATGACCG
63▶ L A W V W V R G L D E L Y A E W S E V V S T N F R D A S G P A M T

7301 **XbaI (7391)** AGATCGGCGAGCAGCCGTGGGGCGGGAGTTCCGCCTGCGCGACCCGGCCGCAACTGCGTGCACCTTTGTGGCAGAGGAGCAGGACTAAATCTAGAATTA
96▶ E I G E Q P W G R E F A L R D P A G N C V H F V A E E Q D •

7401 TCCCTAATACCTGCCACCCACTCTTAATCAGTGGTGAAGAACGGTCTCAGAAGTCTTTGTTTCAATTGGCCATTTAAGTTTAGTAGTAAAAGACTGGT

7501 TAATGATAACAATGCATCGTAAACCTCAGAAGGAAAGGAGAATGTTTTGTGGACCCTTTGGTTTTCTTTTTGCGTGTGGCAGTTTTAAGTTATTAG

7601 **ScaI (7614)** TTTTTAAATCAGTACTTTTTAATGGAAACAATTTGACCAAAAATTTGTACAGAATTTTGGAGCCATTAAAAAGTTAAATGAGAAACCTGTGTGTTCC
7701 CTTTGGTCAACACCGAGACATTTAGGTGAAAGACATCTAATTCTGTTTTACGAATCTGGAAACTTCTTGAATGTAATTCTTGAGTTAACACTTCTGG
7801 GTGGAGAATAGGGTTGTTTTCCCCACATAATTGGAAGGGGAAGGAATATCATTTAAAGCTATGGGAGGTTGCTTTGATTACAACACTGGAGAGAAAT
7901 GCAGCATGTTGCTGATTGCCTGTCACTAAAACAGGCCAAAACCTGAGTCTTGGGTTGCATAGAAAGCTG