

pTRIOZ-hIgG1L2

Plasmid for high yield production of recombinant human IgG1 lambda 2 mAbs

Catalog code: ptrioz-higg1l2

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

For research use only

Version 22B14-MM

PRODUCT INFORMATION

Contents

- 20 µg of pTRIOZ-hIgG1L2 plasmid provided as lyophilized DNA
- 1 ml of Zeocin® (100 µg/ml)

Storage and Stability

- pTRIOZ-hIgG1L2 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-hIgG1L2 expresses the constant region of the heavy (CH) chain from human IgG1 and the constant region of the human lambda 2 light chain (CL). pTRIOZ-hIgG1L2 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, Acc65I, and AvrII* -3'.
- **Human λ2 light chain:** The constant region of the human lambda 2 light chain (IGLC2).
- **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-1alpha 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-hIgG1L2

- **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 Asia: +852 3622-3480
E-mail: info@invivogen.com



PLASMID RESUSPENSION

- Centrifuge the tube containing the lyophilized pTRIOZ-hlgG1L2 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-hlgG1L2, the constant region of the human IgG1 heavy chain is preceded by a MCS containing four unique 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-hlgG1L2, *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

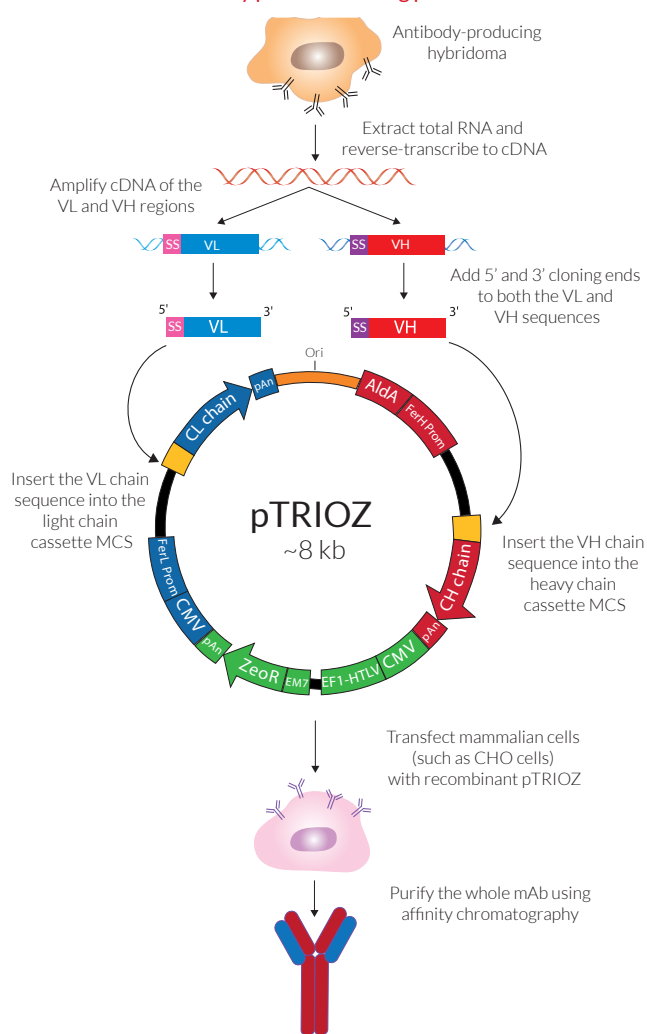
In pTRIOZ-hlgG1L2, the constant region of the human lambda 2 light chain is preceded by a MCS containing six restriction sites: *SgrAI*, *Ascl*, *PmeI*, *NcoI*, *Acc65I* and *AvrII*. We recommend using the *SgrAI* restriction site for insertion of the 5' end of the mAb VL chain (including the native signal sequence).

In pTRIOZ-hlgG1L2, *AvrII* 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 *AvrII* site at the 3' end of the VL chain, in frame with the constant region of the human lambda 2 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.

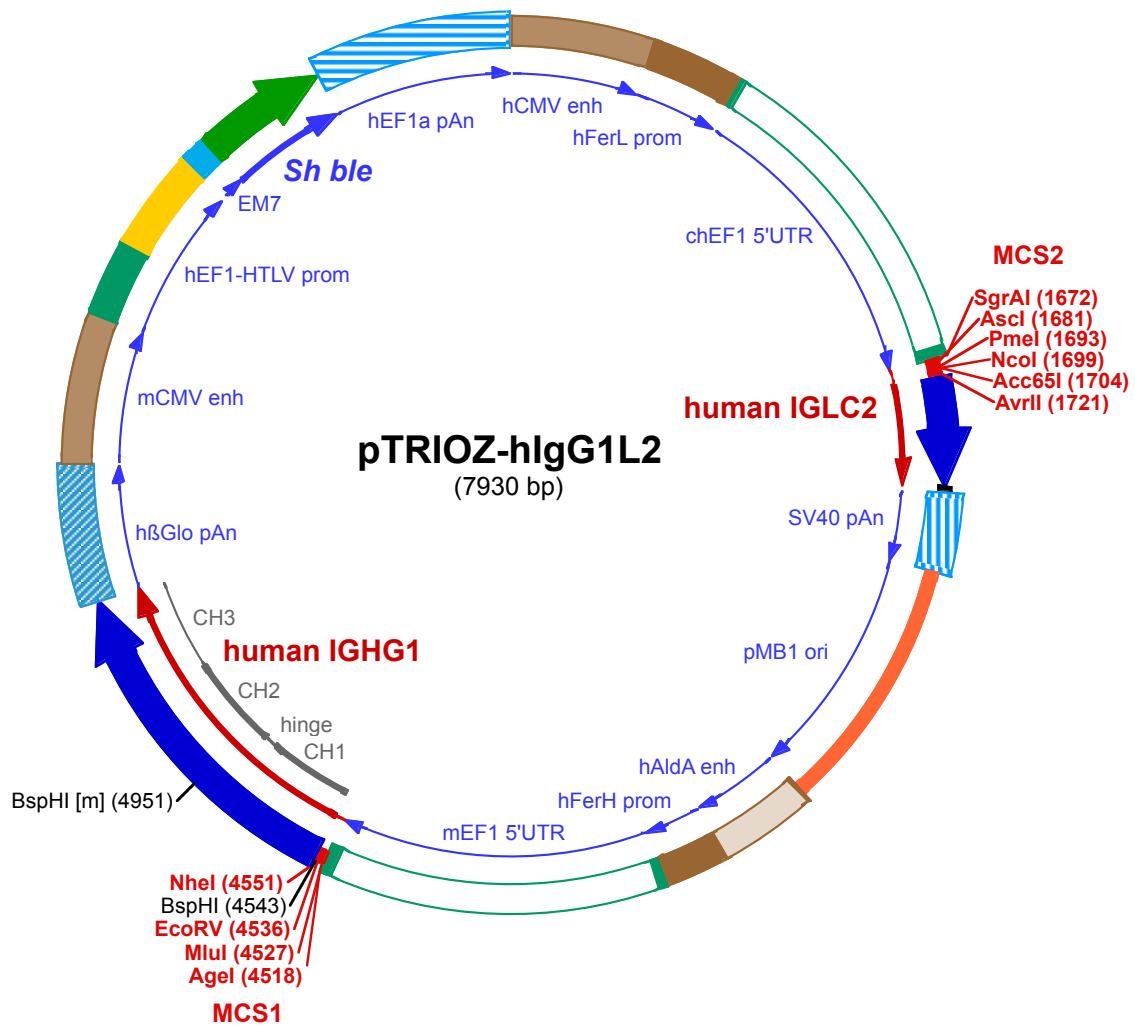
1. Prentice, H.L. et al., 2007. High level expression of proteins using sequences from the ferritin heavy chain gene locus. *J Biotech.* 128:50-60. 2. Rita costa, A. et al., 2010. Guidelines to cell engineering for monoclonal antibody production. *Eur J Pharm Biopharm.* 74(2):127-138.

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 Asia: +852 3622-3480
E-mail: info@invivogen.com



1 CCTGCAGGCGTTACATAACTTACGGTAAATGGCCCGCCTGGCTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAA

101 CGCCAATAGGACTTTCCATTGACGTCAATGGGTGGAGTATTTACGGTAAACTGCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCC

201 TATTGACGTCAATGACGGTAAATGGCCCGCCTGGCATTATGCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATC

301 GCTATTACCATGATGATGCGGTTTTGGCAGTACATCAATGGCGTGGATAGCGGTTTGACTCACGGGGATTCCAAGTCTCCACCCATTGACGTCAATG

401 GGAGTTTGTGTTGACTAGTCAGGGCCCAACCCCCCAAGCCCCATTTCACAACACGCTGGCGCTACAGGCGGTGACTTCCCTTGCTTTGGGGCGGG

501 GGGCTGAGACTCTATGTGCTCCGATTGGTCAGGCACGGCCTTCGGCCCCGCTCCTGCCACCGCAGATTGGCCGCTAGGCCTCCCCGAGCGCCTGCC

601 TCCGAGGGCCGGCGACCATAAAAGAAGCCGCCCTAGCCACGTCCCCTGCAGTTCGGCGGTCCCGGGTCTGTCTCAAGCTTGCCGCGAGAACAAGCAGg

701 taagtgccgtgtgtggttcccgcgggcctggcctctttacgggttatggccttgcgtgccttgaattacttccatgcccctggctgcagtacgtgat

801 ttgatcccagccttcgggttgaagtgggtgggagagttcgaggccttgcgcttaaggagccccttcgctcgtgcttgagttgaggcctggcttggcg

901 ctggggccgcccgtgctaactcgttggcaccttcgcgctgtctcgtgctttcgttaagtctctagccattaaaatTTTTgataaccagctgcgagc

1001 cttttttctgpcgagatagcttctaataatgcccagcagatctgcacactggtatctcggtttttggggccgcccggcgagcggggcccgtgcgtccc

1101 agcgcacatgttcggcgaggcggggcctgcgagcggccaccgagaatcggagcggggtagtctcaactggccggcctgctcgttgcctggcctgcg

1201 gccgcccgtatcggcccctggcggaaggctggcccggctcggcaccagttgcgtgagcggaaagatggccgcttcccggcccctgctgcagggagc

1301 tcaaatggaggacgcccgggagagcgggggggtgagtcaccacacaaagaaaaggccttctcctcatccgtcgttcatgtgactcca

1401 cggagtagcggcgcccgtccaggcacctcgattagttgctgagcttttgagtagctcgtccttaggttgggggaggggtttatgcatggagtttcc

1501 ccacactgagtggggtggagactgaagagttaggccagcttggcacttgatgtaattctccttgaatttggcctttttgagtttggatcttgcctcattc

1601 tcaagcctcagacagtggttcaagttttttcttccatttcagGTGTCGTGAAAAC TACCCCTAAAAGCCACCGGCGAGGCGGCCAAGTTTAAACACC

SgrAI (1672) Ascl (1681) PmeI (1693) NcoI (1699)

1701 ATGGGTACCAAGCTTACCGTCTTAGGTCAGCCCAAGGCTGCCCTCGGTCACTCTGTCCCGCCTCCTCTGAGGAGCTTCAAGCCAACAAGCCACAC

1▶ G T K L T V L G Q P K A A P S V T L F P P S S E E L Q A N K A T

1801 TGGTGTGTCTATAAGTACTTCTACCCGGGAGCCGTGACAGTGGCCTGGAAGGCAGATAGCAGCCCCGTCAAGGCGGGAGTGAGACCACCACCCCTC

26▶ L V C L I S D F Y P G A V T V A W K A D S S P V K A G V E T T T P S

1901 CAAACAAAGCAACAAGTACGCGCCAGCAGCTATCTGAGCCTGACGCCTGAGCAGTGAAGTCCCACAGAAGTACAGCTGCCAGGTCACGCATGAA

59▶ K Q S N N K Y A A S S Y L S L T P E Q W K S H R S Y S C Q V T H E

2001 GGGAGACCCTGGAGAAGACAGTGGCCCTACAGAATGTTTCATAGCTAGGACACTAGTCACCTAGCTGGCCAGACATGATAAGATAACATTGATGAGTTG

93▶ G S T V E K T V A P T E C S •

2101 GACAAACCACAAC TAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGT

2201 TAACAACAACAATTGCATTCATTTTATGTTTCAGGTTCAAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATG

2301 TTAATTAAGTACCATGACCAAAATCCCTTAACGTGAGTTTTTCGTTCCACTGAGCGTCAGACCCCGTAGAAAAGATCAAAGGATCTTCTTGAGATCCTTT

2401 TTTTCTGCGGTAATCTGCTGCTTGCAACAAAAAACCACCGCTACCAGCGTGGTTTGTGTTGCCGATCAAGAGCTACCAACTCTTTTTCCGAAGGTA

2501 ACTGGCTTCAGCAGAGCGCAGATACCAAACTGTTCTTCTAGTGTAGCCGTAGTTAGGCCACCCTCAAGAAGTCTGTAGCACCCTACATACTCTG

2601 CTCTGCTAATCCTGTTACCAGTGGCTGCTGCCAGTGGCGATAAGTCTGTCTTACCGGTTGGACTCAAGACGATAGTTACCAGATAAGGCGCAGCGGTC

2701 GGGCTGAACGGGGGTTCTGTGCACACAGCCAGCTTGGAGCGAACGACCTACCCGAAGTACGATACCTACAGCGTGAAGTATGAGAAAAGCGCCACGCTT

2801 CCCGAAGGGAGAAAAGCGGACAGGTATCCGGTAAGCGGAGGGTCCGAACAGGAGAGCGCACGAGGGAGCTTCCAGGGGAAACGCCTGGTATCTTTATA

2901 GTCCTGTGCGGTTTCCGCACCTGACTTGGAGCGTCAATTTTGTGATGCTGTCAGGGGGCGGAGCCTATGAAAAACGCCAGCAACGCGGCTTTTT

3001 ACGGTTCTGGCCTTTTGTGCGCTTTTGTCCATGTTCTTAATTAACCTGCaggcgaactcagcttctccttgcgtttccgacttttccatccgctcct

3101 ccacttccccgttccgcctccccattgccaacattctggctgagtcacggcgccccagagcgcgccaggctgggggaaaggagcagaaggaggggccc
3201 tagcgaccgcgggatgtggtccgagtcacgtccgaggggggtggggaggatcggttctcggcgccccttctagcgcggcctctgggtcgcg
3301 ctctcggggcgcccgtagcccagtcctgctgACTAGTTCGCCAGAGCGCGGAGGGCTCCAGCGGCCCCCTCCCCACAGCAGGGCGGGTCCC
3401 GCGCICACCGGAAGGAGCGGGCTCGGGGCGGGCGCGCTGATTGGCCGGGGCGGGCTGACGCCGACGCGGTATAAGAGACCACAAGCGACCCGAGGG
3501 CCAGACGTTCTTCGCCAAGCTTGCCGTGAGAACGACGgtgagggcggggtggttccgcgggcccagctggaggctctgctcagagcgggccc
3601 gccccgtgtcgtcggcggggatttagctgcgagcattcccgttcgagttcggggcgcgggaggcagagtgcgaggcctagcggcaacccgtagcc
3701 tcgctcgtgtccggcttgaggcctagcgtggtgtccgcgccgcccgcgtgctactccggcgcactctggtctttttttttttgttgttgttggcc
3801 tgctgccttcgattgccgttcagcaataggggtaacaaggagggtgctggggcttctcgcggagcccggagaggtcatggttggggaggaatgga
3901 gggacaggagtggcggctggggcccgccttcggagcacatgtccgacgccacctggatggggcaggcctggggttttccgaagcaaccaggct
4001 ggggttagcgtgccgaggccatgtggccccagacccggcacgatctggcttggcggcgccgcttgcctcctcctaactagggtgaggccatcccg
4101 tccggcaccagttgcgtcgtggaagatggcgcctccgggcccctgttgcaaggagctcaaagtggaggacgcggcagcccgggtggagcgggctga
4201 gtacccacacaaaggaagggcctggtccctaccggctgctgcttctgtgaccccggtgctctatcggccgcaatagtcacctcgggcttttagc
4301 acggctagtcgcggcgggggaggggatgtaatggcgttggagtttggcatttgggtgggagactagttagccagcctggcgtggaagtcat
4401 tttggaatttgtccccttgagttttgagcggagctaattctcgggcttcttagcggttcaaaggatcttttaaaccttttttaggTGTTGTGAAAAC

4501 ACCGCTAATTCAAAGCAACCGGTCGCACGCGTAGATATCACGTCATGAAAGCTAGCACCAAGGGCCATCGGTCTTCCCCCTGGCACCTCCTCCAAGAG
A S T K G P S V F P L A P S S K S

4601 CACCTCTGGGGCACAGCGCCCTGGGCTGCCTGGTCAAGGACTACTTCCCCGAGCCGGTACGGTGTCTGGAAGTCAAGGCGCCCTGACCAGCGGGCTG
17> T S G G T A A L 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

4701 CACACCTTCCCGCTGTCTACAGTCTCAGGACTCTACTCCCTCAGCAGCGTGGTACCGTGCCTCCAGCAGCTTGGGCACCCAGACCTACATCTGCA
51> H T F P A V L Q 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

4801 ACGTGAATCACAAGCCAGCAACACCAAGTGGACAAGAAAGTTGAGCCAAATCTTGTGACAAAAGTACACATGCCACCGTGCCAGCACCTGAACT
84> N V N H K P S N T 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

4901 CTGGGGGACCCTCAGTCTTCTTCCCCCAAACCAAGGACACCTCATGATCTCCGGACCCCTGAGGTCACATGCGTGGTGGTGGACGTGAGC
117> L G G P S V F L 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

5001 CACGAAGACCCTGAGGTCAAGTCAACTGGTACGTGGACGGCTGGAGGTGCATAATGCCAAGACAAAGCCGCGGGAGGAGCAGTACAACAGCACGTACC
151> H E D P E V K F 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

5101 GTGTGGTACAGCTCTCACCGTCTGCACCAGGACTGGCTGAATGGCAAGGAGTACAAGTGAAGGTCTCAAACAAAGCCCTCCAGCCCCATCGAGAA
184> R V V S V L T V L 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

5201 AACCATCTCAAAGCCAAAGGGCAGCCCGAGAACCACAGGTGTACACCCTGCCCCATCCCGGGACGAGCTGACCAAGAACCAGGTGAGCTGACCTGC
217> T I S K A K G Q 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

5301 CTGGTCAAAGGCTTCTATCCCAGCGACATCGCGTGGAGTGGGAGAGCAATGGCAGCCGGAGAACAACACTACAAGACCAGCCTCCCGTGTGGACTCCG
251> L V K G F Y P S 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

5401 ACGGCTCTTCTTCTCTACAGCAAGCTCACCGTGGACAAGAGCAGGTGGCAGCAGGGGAACGCTTCTCATGCTCCGTGATGCATGAGGCTCTGCACAA
284> D G S F F L Y S K 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

5501 CCACTACAGCAGAAGAGCCTCTCCCTGTCTCCGGTAAATGAGTCTAGAAGCTCGCTTTCTTGCTCCAATTTCTATTAAGGTTCTTTGTTCCCT
317> H Y T Q K S L S L S P G K •

5601 AAGTCCAACACTAAACTGGGGATATTATGAAGGGCCTTGGAGCATCTGGATTCTGCCTAATAAAAAACATTTATTTTCATTGCAATGATGATTTAAAT
5701 TATTTCTGAATATTTTACTAAAAAGGAATGTGGGAGTCAAGTTCATTTAAAAACATAAAGAAATGAAGAGCTAGTTCAAACCTTGGGAAAATACACTATA

5801 TCTTAACTCCATGAAAGAAGGTGAGGCTGCAACAGCTAATGCACATTGGCAACAGCCCTGATGCCTATGCCTTATTCATCCCTCAGAAAAGGATTCA

5901 AGTAGAGGCTTGATTTGGAGGTTAAAGTTTTGCCATGCTGTATTTTAGAATTCTGCAGGAGTCAATGGGAAAAACCCATTGGAGCCAAGTACACTGACT
6001 CAATAGGGACTTTCCATTGGGTTTTGCCAGTACATAAGGTCAATAGGGGGTGAAGTCAACAGGAAAGTCCCATTGGAGCCAAGTACATTGAGTCAATAGG
6101 GACTTTCCAATGGGTTTTGCCAGTACATAAGGTCAATGGGAGGTAAGCCAATGGGTTTTCCATTACTGCAIGTATACTGAGTCATTAGGGACTTTC
6201 CAATGGGTTTTGCCAGTACATAAGGTCAATAGGGGTGAATCAACAGGAAAGTCCCATTGGAGCCAAGTACACTGAGTCAATAGGGACTTTCCATTGGGT
6301 TTTGCCAGTACAAAAGGTCAATAGGGGGTGAAGTCAATAGGGTTTTCCATTATTGGCACATACATAAGGTCAATAGGGGTGACTAGTCAGTGGGCAGAG
6401 CGCACATCGCCCGGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTAGAGAAGGTGGCGCGGGTAAACTGGGAAAGTGATGTCGTGACTG
6501 GCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCCGTGAACGTTCTTTTTCGAACGGGTTTCCGCCAGAACACAGCTG
6601 AAGCTTCGAGGGCTCGCATCTCTCTTACGCGCCGCCCTACCTGAGGCCCATCCACGCCGTTGAGTCCGCTTCTGCCGCTCCCGCTGTG
6701 GTGCCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGTCAAGTGCAGACCGGGCCTTTGTCCGGCGCTCCCTTGGAGCTACCTAGACTCAGCCG
6801 GCTCTCCACGCTTTCCTGACCCTGCTGCTCAACTCTACGCTTTTGTTCGTTTTCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGCGCTACA
6901 AACAGTAGTTGACAATTAATCATCGGCATAGTATATCGGCATAGTATAATACGACTCACTATAGGAGGGCCATCATGGCCAAGTTGACCAGTCCGTTCC
7001 GGTGTCACCGCGCGAGCTCGCCGGAGCGGTGAGTTCTGGACCGACCGGCTCGGGTTCTCCCGGACTTCGTGGAGGACGACTTCGCTGGTGTGGTC
9▶ V L T A R D V A G A V E F W T D R L G F S R D F V E D D F A G V V
7101 CGGACGACGTGACCCTGTTTCATCAGCGCGGTCCAGGACCAGGTGGTCCCGGACAACACCCTGGCCTGGGTGTGGGTGCGCGGCTGGACGAGCTGTACG
43▶ R D D V T L F I S A V Q D Q V V P D N T L A W V W V R G L D E L Y
7201 CCGAGTGGTGGAGGTGCTGTCACGAACCTCCGGGACGCCTCCGGGCGGCCATGACCGAGATCGGCGAGCAGCCGTGGGGCGGGAGTTCGCCCTGCG
76▶ A E W S E V V S T N F R D A S G P A M T E I G E Q P W G R E F A L R
7301 CGACCGGCGGCAACTGCGTCACTTTGTGGCAGAGGAGCAGGACTAACTAGAATTATCCCTAATACCTGCCACCCCACTCTTAATCAGTGGTGAA
109▶ D P A G N C V H F V A E E Q D •
7401 GAACGGTCTCAGAAGTGTGTTTCAATTGGCCATTTAAGTTTAGTAGTAAAAGACTGGTTAATGATAACAATGCATCGTAAAACCTCAGAAGGAAAGG
7501 AGAATGTTTTGTGGACCACTTTGGTTTTCTTTTTGCGTGTGGCAGTTTTAAGTTATTAGTTTTAAAATCAGTACTTTTTAATGAAACAACCTTGACCA
7601 AAAATTTGTCACAGAATTTTGGACCCATTAATAAAGTTAAATGAGAAACCTGTGTTCCTTTGGTCAACACCGAGACATTTAGGTGAAAGACATCTAA
7701 TTCTGTTTTACGAATCTGAAACTTCTGAAATGTAATTCTTGAGTTAACACTTCTGGGTGGAGAATAGGGTTGTTTTCCCCCACATAATTGGAAGG
7801 GGAAGGAATATCATTTAAAGCTATGGGAGGTTGCTTTGATTACAACACTGGAGAGAAATGCAGCATGTTGCTGATTGCTCACTAAAACAGGCCAAA
7901 AACTGAGTCCTTGGGTTGCATAGAAAGCTG
▶