

pFUSEss-CHlg-hD

Plasmid featuring the constant region of the human IgD (allele 2) heavy chain, and the IL2 signal sequence

Catalog code: pfusess-hchd2

For research use only

Version 24J28-MM-v37

PRODUCT INFORMATION

Content:

- 20 µg of pFUSEss-CHlg-hD plasmid provided as lyophilized DNA
- 1 ml of Zeocin® (100 µg/ml)

Storage and Stability:

- Product is shipped at room temperature.
- Lyophilized DNA should be stored at -20°C and is stable 3 months.
- Resuspended DNA should be stored at -20°C and is stable up to 1 year.
- Store Zeocin® at 4 °C or at -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.

Materials required for antibody generation & isotype switching

- pFUSE2ss-CLlg plasmid that features the constant region of the kappa or lambda light chains. pFUSE2ss-CLlg plasmids are selectable with blasticidin (sold separately, see RELATED PRODUCTS).
- pFUSEss-CHlg plasmid for the constant region of the heavy chain. This plasmid is selectable with Zeocin®.

GENERAL PRODUCT USE

pFUSEss-CHlg and pFUSE2ss-CLlg plasmids are designed to change a monoclonal antibody from one isotype to another, therefore, enabling the generation of antibodies with the same antigen affinity but with different effector functions (increased or reduced ADCC and CDC). Furthermore, they can be used to produce entire antibodies from Fab or scFv fragments.

pFUSEss-CHlg and pFUSE2ss-CLlg express the constant regions of the heavy (CH) and light (CL) chains, respectively. They contain a multiple cloning site (MCS) upstream of these constant regions to enable the cloning of the variable (VH and VL) regions of a given antibody. Transfection of mammalian cell lines with the recombinant pFUSEss-CHlg and pFUSE2ss-CLlg pair allows to generate an antibody that can be purified from the supernatant using the appropriate affinity chromatography.

Features of pFUSEss-CHlg and pFUSE2ss-CLlg plasmids

- **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1α (EF-1α) core promoter¹ and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat². 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.
- **MCS:** The multiple cloning site contains several restriction sites that are compatible with many other enzymes, thus facilitating cloning.
- **SV40 pAn:** the Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA³.
- **ori:** a minimal *E. coli* origin of replication to limit vector size, but with the same activity as the longer Ori.
- **CMV enh / hFerL prom:** This composite promoter combines the human cytomegalovirus immediate-early gene 1 enhancer and the core promoter of the human ferritin light chain gene. This ubiquitous promoter drives the expression of the Zeocin®-resistance gene in mammalian cells.
- **IL2 ss:** The human IL2 signal sequence contains 20 amino acids (MYRMQLLSICIALSLALVTNS) and share common characteristics with signal peptides of other secretory proteins. The intracellular cleavage of the IL2 signal peptide occurs after Ser20 and leads to the secretion of the immunoglobulin chain.
- **EM2KC** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*. EM2KC is located within an intron and is spliced out in mammalian cells.
- **βGlo pAn:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription⁴.

pFUSEss-CHlg-hD specific features

- **Human IGHD (IgD allele 2 heavy chain constant region):** When cloning your heavy chain variable region of choice in the MCS, care must be taken to insert the gene in-frame and to preserve the integrity of the heavy chain constant region.
- **Zeo:** Resistance to Zeocin® is conferred by the *Sh ble* gene from *Streptoalloteichus hindustanus*. The same resistance gene confers selection in both mammalian cells and *E. coli*.

References:

1. Kim DW, *et al.* 1990. Use of the human elongation factor 1 alpha promoter as a versatile and efficient expression system. *Mol Cell Biol.* 9(2):217-23.
2. Takebe Y, *et al.* 1988. SR alpha promoter: an efficient and versatile mammalian cDNA expression system composed of the simian virus 40 early promoter and the R-U5 segment of human T-cell leukemia virus type 1 long terminal repeat. *Mol Cell Biol.* 8(1):466-72.
3. Carswell S, & Alwine JC. 1989. Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. *Mol Cell Biol.* 9(10):4248-58.
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 Asia: +852 3622-3480
E-mail: info@invivogen.com



PROTOCOL

Obtaining VH and VL sequences

To obtain the cDNA sequence of the VH and 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. Alternatively 5' RACE can be used. The resulting amplicons must be sequenced.

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.

Cloning into pFUSEss-CHIg and pFUSE2ss-CLIg

Once the VH and VL sequence are known, inserts for cloning into the plasmids can be generated. In pFUSEss-CHIg-hD, the constant region of the human IgD heavy chain is preceded by a multiple cloning site containing five unique restriction sites: EcoRI, EcoRV, XhoI, NheI, and Eco47III. Using EcoRI as the 5' cloning site ensures that the cloned VH will follow the hIL2 signal sequence without unwanted additional amino-acids. Use Eco47III as the 3' cloning site in order to preserve the exact IgD amino acid sequence. We recommend introducing by PCR the EcoRI and Eco47III sites at the VH boundaries. Care must be taken to preserve the correct reading frame when placing the cloning sites on the variable sequence.

Note:

When generating the insert for VL, a *BsiWI* (*pFUSE2ss-CLIg-hk*; human kappa), or *AvrII* (*pFUSE2ss-CLIg-hl2*; human lambda 2) site must be introduced at the 3' end.

Antibody production

Cotransfect mammalian cells, such as 293 and CHO cells, with the recombinant plasmids pFUSE2ss-CLIg encoding the light chain and pFUSEss-CHIg encoding the heavy chain. Antibody production depends greatly on the ratio of heavy chain and light chain expression. Typically, pFUSEss-CHIg to pFUSE2ss-CLIg ratio of 2:3 is used to cotransfect mammalian cells. Since both plasmids share the same plasmid backbone, the appropriate heavy chain to light chain ratio can be easily determined by varying the quantities of plasmids.

OR

Transfect cells using a transfection agent, such as LyoVec™, with the plasmid coding for light chain and select the best clone. Following selection of the best clone, the plasmid coding for the heavy chain clone can be transfected into this clone.

Use blasticidin and Zeocin® to select pFUSE2ss-CLIg and pFUSEss-CHIg respectively.

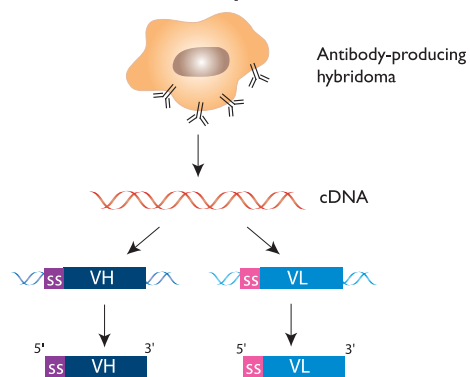
Antibody production can be analyzed by different techniques including SDS-PAGE, flow cytometry, ELISA, or a bioactivity assay.

Antibody purification

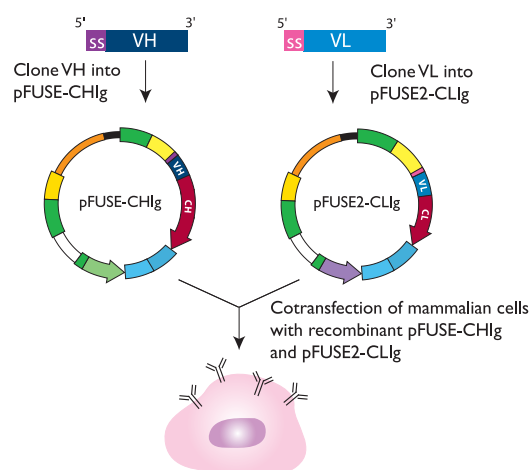
The resulting IgG antibody can be purified from the supernatant using the appropriate Protein A, Protein G or Protein L affinity chromatography.

Antibody generation using pFUSE-CHIg & pFUSE-CLIg

1- Obtention of VH and VL sequences



2- Cloning into pFUSE-CHIg and pFUSE-CLIg



RELATED PRODUCTS

| Product | Catalog Code |
|---------------------|----------------|
| pFUSE2ss-CLIg-hk | pfuse2ss-hclk |
| pFUSE2ss-CLIg-hl2 | pfuse2ss-hcll2 |
| LyoVec™ | lyec-12 |
| Protein L / Agarose | gel-protl-2 |
| Zeocin® | ant-zn-1 |

TECHNICAL SUPPORT

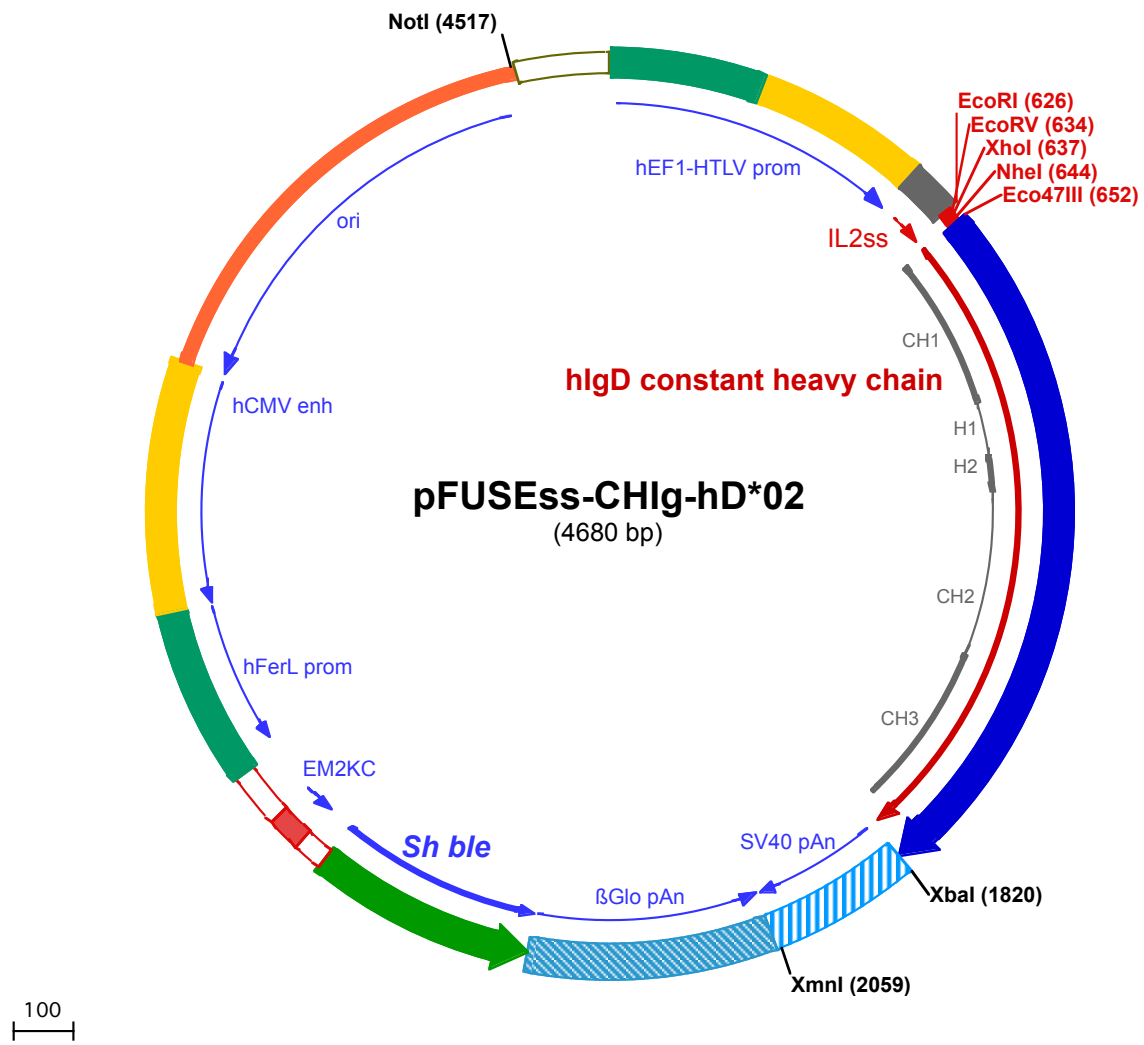
InvivoGen USA (Toll-Free): 888-457-5873

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 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
101 GAGAAGGTGGCGCGGGGTAAACTGGGAAAGTGATGCTGTACTGGCTCCGCCTTTTTCCCGAGGGTGGGGGAGAACCGTATATAAGTGCAGTAGTCGCC
201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCCGCCCTACCTGAGGCC
301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCTGACCTGCTTGTCTAACTCTACGCTTTTGTTCGTTT
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGCGAAGGAGGGCCACCATGTACAGGATGCAACTCCTGTCTTGA
1 M Y R M Q L L S C
EcoRV (634) NheI (644)
EcoRI (626) XhoI (637) Eco47III (652)
601 TTGCACTAAGTCTTGCACTTGTACGCAATTGATATCTCGAGTCTAGCAGCGCTCCACCAAGGCTCCGGATGTGTTCCCATCATATCAGGGTGCAGA
10 I A L S L A L V T N S 1 A P T K A P D V F P I I S G C R
701 CACCCAAAGGATAACAGCCCTGTGGTCTGGCATGCTTGATAACTGGGTACCACCAACGCTCCGTGACTGTACCTGGTACATGGGGACAGAGCCAGC
17 H P K D N S P V V L A C L I T G Y H P T S V T V T W Y M G T Q S Q
801 CCCAGAGAACCTCCCTGAGATACAAAGACGGGACAGCTACTACATGACAAGCAGCCAGCTCTCCACCCCTCCAGCAGTGGCGCAAGGCGAGTACAA
50 P Q R T F P E I Q R R D S Y Y M T S S Q L S T P L Q Q W R Q G E Y K
901 ATGCGTGGTCCAGCACACCGCCAGCAAGAGTAAGAAGGAGATCTTCCGCTGGCAGAGTCTCAAAGGCACAGCCTCCTCAGTCCCCTGCACAACCC
83 C V V Q H T A S K S K K E I F R W P E S P K A Q A S S V P T A Q P
1001 CAAGCAGAGGGCAGCCTGCCAAGGCAACCACAGCCCCAGCCACCACCGTAACACAGGAAGAGGAGGAGAAGAGAAGAAGGAGAAGGAGAAAGAGG
117 Q A E G S L A K A T T A P A T T R N T G R G G E E K K K E K E K E
1101 AACAGAAGAGAGAGACAAAGACACCAGAGTGTCCGAGCCACCCAGCCTTTGGCGTCTACCTGCTAACCCCTGCAGTGCAGGACCTGTGGCTCCG
150 E Q E E R E T K T P E C P S H T Q P L G V Y L L T P A V Q D L W L R
1201 GGACAAAGCCACCTTACCTGCTTCTGGTGGGCGAGTACCTGAAGGATGCTACCTGACCTGGGAGGTGGCTGGGAAGTCCCCACAGGGGGCGTGGAG
183 D K A T F T C F V V G S D L K D A H L T W E V A G K V P T G G V E
1301 GAAGGGCTGTGGAGCGGCACAGCAACGGTCCCAGAGCCAGCACAGCCGTGACCCTGCCAGGTCTTGTGGAACCGGGGACCTCCGTACCTGCA
217 E G L L E R H S N G S Q S Q H S R L T L P R S L W N A G T S V T C
1401 CACTGAACCATCCAGCCTCCACCCAGAGGTTGATGGCGTGGAGAACCCGCTGCGCAGGCACCCGTCAAGTCTTCCCTGAACCTGTGGCTCGTC
250 T L N H P S L P P Q R L M A L R E P A A Q A P V K L S L N L L A S S
1501 TGACCCTCCCAGGGCGCCTCGTGGCTCCTGTGTGAGGTGTCTGGCTTCTGCCCCCAACATCCTCCTGATGTGGCTGGAGGACCAGCGTGAAGTGAAC
283 D P P E A A S W L L C E V S G F S P P N I L L M W L E D Q R E V N
1601 ACTTCTGGGTTTGCCCCGCACGCCCCCTCCACAGCCAGGAGCACCAGTTCTGGGCTGGAGTGTGCTGCGTGTCCAGCCCCGCCAGCCCTCAGC
317 T S G F A P A R P P P Q P R S T T F W A W S V L R V P A P P S P Q
1701 CAGCCACCTACACGTGTGGTTCAGCCACGAGGACTCCCGACTCTGCTCAACGCCAGCCGGAGCCTAGAAGTCAAGTATGTAACAGACCATGGCCCCAT
350 P A T Y T C V V S H E D S R T L L N A S R S L E V S Y V T D H G P M
XbaI (1820)
1801 GAAATGATCCCGACCAGATCTAGACCTAGTGGCCAGACATGATAAGATACATTGATGAGTTTGACAAACCACAACCTAGAATGCAGTGAAAAAATGC
383 K •
1901 TTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACATTCATTCTTTATGTTTCAGG
XmnI (2059)
2001 TTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTAATTCTAAAATACAGCATAGCAAACTTTAACCTCC
2101 AAATCAAGCCTCTACTGAACTCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTGCAATGTGCATTAGCTGTTTGCAGCCTCACCTCT
2201 TTCATGGAGTTTAAAGATATAGTGTATTTTCCCAAGGTTTGAAGTACTCTTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTAGT
2301 AAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAG
2401 TTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCTCAGTCTGCTCTGCCACAA
125 D Q E E A V F
2501 AGTGCACGAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCCCGCCCCACGGCTGCTCGCGATCTCGGTTCATGGCCGGCCGGAGGCGTCCCGGAAGTT
117 H V C N G A P D R L A F E R G W P Q E G I E T M A P G S A D R F N

2601 CGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCCGCGCACCCACCCAGGCCAGGGTGTGTCCGGCACCACCTGGTCCTGGACCGCG
84 T S V V E S W E A Y L E D L G R V W V W A L T N D P V V Q D Q V A
2701 CTGATGAACAGGGTCACGTCGTCCCGGACCACACCGCGAAGTCGTCTCCACGAAGTCCCGGAGAACCCGAGCCGGTCCGGTCCAGAACTCGACCGCTC
50 S I F L T V D D R V V G A F D D E V F D R S F G L R D T W F E V A G
2801 CGGCGACGTCGCGCGGGTGAAGCACCAGGACGGCACTGGTCAACTTGGCCATGATGGCTCTCctgtcaggagaggaagagaagaaggttagtacaatt
17 A V D R A T L V P V A S T L K A M
2901 gCTATAGTGAGTTGATTATACTATGCAGATATACTATGCCAATGATTAATTGCAAAGTGGCTGCAgggttcatagtgccacttttcctgcactgcc
←
3001 ccatctcctgccaccctttccaggcatagacagtcagtgacttacCAAACACAGGAGGGAGAAGGCAGAAGCTTGAGACAGACCCGCGGGACCGCC
←
3101 GAACTGCGAGGGGACGTGGCTAGGGCGGCTTCTTTTATGGTGCGCCGCCCTCGGAGGCAGGGCGCTCGGGGAGGCCCTAGCGGCAATCTGCGGTGGCAG
←
3201 GAGGCGGGGCCGAAGGCCGTGCCTGACCAATCCGGAGCACATAGGAGTCTCAGCCCCCGCCCAAAGCAAGGGGAAGTCACGCGCTGTAGCGCCAGCG
←
3301 TGTTGTGAAATGGGGCTTGGGGGGTTGGGGCCCTGACTAGTCAAAACAACTCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCA
←
3401 AACCGTATCCACGCCATTGATGTAAGTCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATA
←
3501 AGGTCATGTAAGTGGCATAATGCCAGGCGGGCATTACCCTGATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGTACTGCAAGTG
←
3601 GGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTACTATGGAAACATACGTCAATTATTGACGTCAATGGGCGGGG
←
3701 TCGTTGGGCGGTGAGCCAGGCGGGCCATTTACCCTAAGTTATGTAACGCCCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAA
←
3801 CCGTAAAAAGGCCGCTTGTGCGTTTTTCCATAGGCTCCGCCCTGACGAGCATCAGAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACA
←
3901 GGACTATAAGATACCAGGCTTTCCCTGGAAGCTCCCTCGTGGCTCTCTGTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTT
←
4001 CGGGAAGCGTGGCGCTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCA
←
4101 GCCCGACCGCTGCGCTTATCCGTAATATCGTCTTGTAGTCCAAACCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGC
←
4201 AGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGC
←
4301 CAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGAAGCAGCAGATTACGCGCAG
←
4401 AAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAA
←
NotI (4517)
4501 TTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTATTACATCTGTGTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCA
4601 AAACAAAACGAAACAAAACAACTAGCAAAATAGGCTGTCCCGAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA