

pFUSE-SEAP-mG1-Fc

Plasmid designed for the expression of a SEAP-Fc Fusion protein

Catalog code: pfuse-mg1sp

<https://www.invivogen.com/pfuse-seap-mg1-fc>

For research use only

Version 22H30-MM

PRODUCT INFORMATION

Contents

- 20 µg of pFUSE-SEAP-mG1-Fc plasmid provided as lyophilized DNA
- 1 ml of Zeocin® (100 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 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 and lyophilized.
- Expression of SEAP-mG3Fc was confirmed by using QUANTI-Blue™ Solution.
- SEAP-mG3Fc protein was purified using protein G affinity chromatography following manufacturer's protocol.

GENERAL PRODUCT USE

pFUSE-SEAP-Fc plasmids express a SEAP-Fc fusion protein generated by fusing the gene encoding for human secreted alkaline phosphatase (SEAP) and the Fc region of an immunoglobulin G (IgG).

pFUSE-SEAP-Fc plasmids yield high levels of Fc-Fusion proteins. The level of expression is usually in the µg/mL range. They can be transfected in a variety of mammalian cells, including myeloma cell lines, Chinese hamster ovary (CHO) cells, monkey COS cells and human embryonic kidney (HEK)293 cells. These cells are commonly used in protein purification systems.

SEAP-Fc fusion proteins are secreted and can be easily detected in the supernatant of pFUSE-SEAP-Fc-transfected cells by using QUANTI-Blue™ Solution, a SEAP detection medium.

WARNING: mIgG1-Fc has a very low affinity to Protein A and Protein G making it difficult to purify.

PLASMID FEATURES

• **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.

• **SEAP-mIgG1-Fc** was generated by fusing the gene encoding for human SEAP with the Fc region of mouse IgG3. This region comprises the CH2 and CH3 domains of the IgG heavy chain and the hinge region. The hinge serves as a flexible spacer between the SEAP and Fc moieties, allowing each part of the molecule to function independently.

• **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.

• **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.

• **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*.

• **βGlo pAn:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription⁴.

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.* 10(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

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 other commonly used laboratory *E. coli* strains, such as DH5α.

Zeocin® usage

This antibiotic can be used for *E. coli* at 25 µg/ml in liquid or solid media and at 50-200 µg/ml to select Zeocin®-resistant mammalian cells.

RELATED PRODUCTS

Product	Catalog Code
LyoVec™	lyec-12
QUANTI-Blue™ Solution	rep-qbs

TECHNICAL SUPPORT

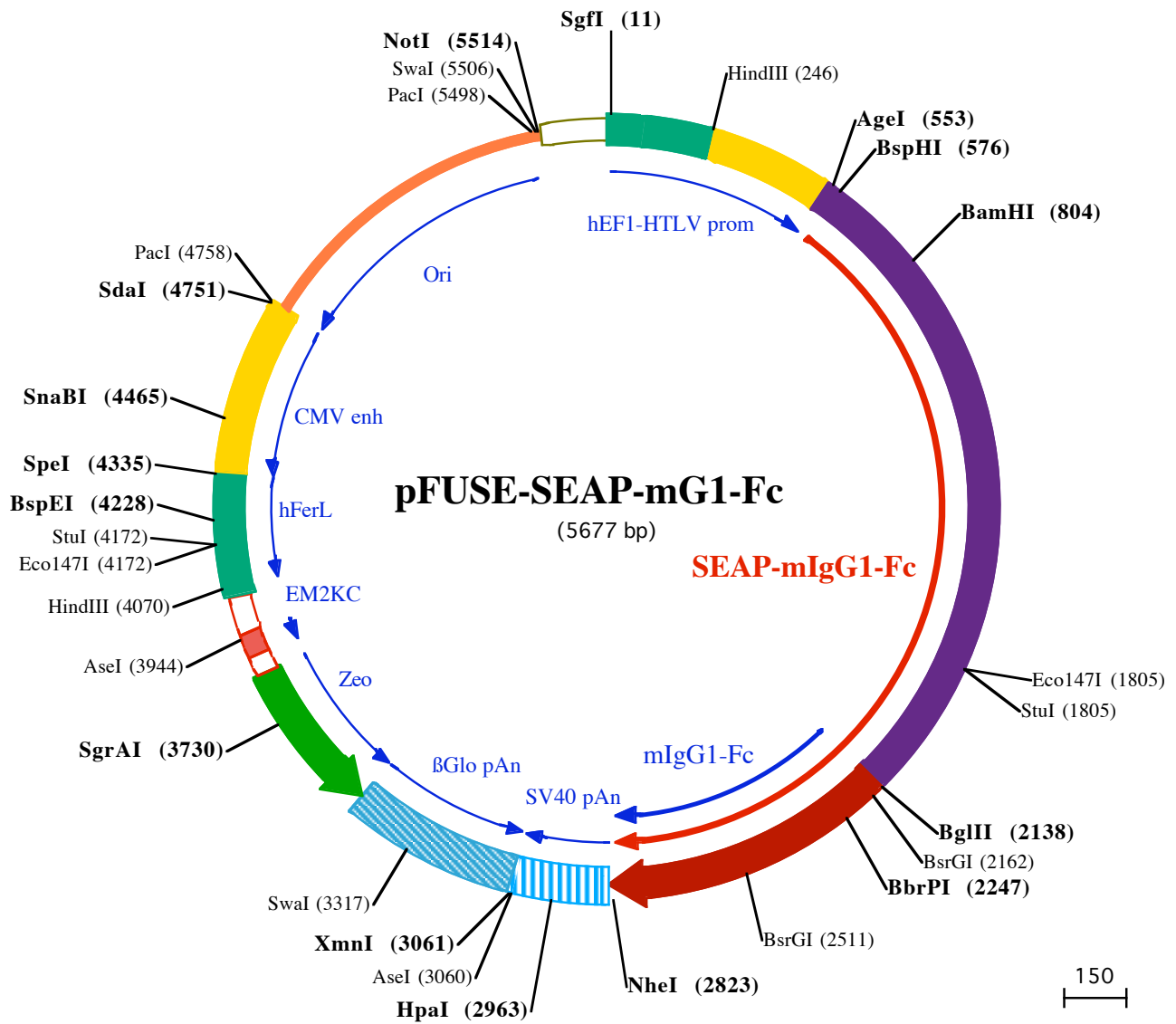
InvivoGen USA (Toll-Free): 888-457-5873

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



SgfI (11)
1 GGATCTGCGATCGCTCCGGTCCCGTCCAGTGGGCGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGTTCGCAATTGAACGGGTGCCTA
101 GAGAAGGTGGCGGGGTAAGTGGGAAAGTGATGTCGTGTAAGTGGCTCCGCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (246)
201 GTGAACGTTCTTTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACCGCGCCCGCCCTACCTGAGGCC
301 GCCATCCACGCGGTTGAGTCGCGTCTGCCGCTCCCGCTGTGGTGCCTCCTGAAGTGCCTCCGCTTAGTAAAGTTAAAGCTCAGGTCGAGACC
401 GGGCCTTTGTCCGCGCTCCCTTGAGGCTACCTAGACTACGCGGCTCTCCAGCTTTGCTGACCTGCTTGTCTCAACTCTACGCTTTGTTTCTGTTT

AgeI (553) BspHI (576)
501 TCTGTTCTGGCGCTTACAGATCCAAGCTGTGACCGGCCCTACCTGAGATCACCGTTTCAGCTGAGGAGGCACATCATGATTCTGGGGCCCTGCATGCT
601 GCTGCTGCTGCTGCTGGCCTGAGGCTACAGTCTCCCTGGGCATCATCCAGTGTAGGAGGAGAACCAGGACTTCTGGAACCGCGAGGCGAGCCGAG
801 CCAAGTCTTAAAGGGCAGAAGAAGGACAACTGGGGCTGAGATACCCCTGGCTATGGACCGCTTCCCATATGTGGCTGTCCAAGACATACAATGT
75▶IaArgIleLeuLysGlyGlnLysLysAspLysLeuGlyProGluIleProLeuAlaMetAspArgPheProTyrValAlaLeuSerLysThrTyrAsnVal
901 AGACAAATATGTCCAGACAGTGGAGCCACAGCCACGCGCTACCTGTGCGGGTCAAGGGCACTTCCAGACCATTTGGCTTGTGAGTGCAGCGCCCGCTTT
108▶IAspLysHisValProAspSerGlyAlaThrAlaThrAlaTyrLeuLysGlyValLysGlyAsnPheGlnThrIleGlyLeuSerAlaAlaAlaArgPhe
1001 AACAGTGAACACGACAGCGCGCAACGAGGTCTCTCGTGTGAATCGGCAAGAAAGCAGGGAAGTCACTGGGAGTGGTAACCACACACGAGTGC
142▶AsnGlnCysAsnThrThrArgGlyAsnGluValIleSerValMetAsnArgAlaLysLysAlaGlyLysSerValGlyValValThrThrArgValG
1101 AGCACGCTCGCCAGCCACCTACGCCACAGGTAACCGCACTGACTCGGACCGCCGCTTTCGAGCTGGAGCCGACGTGCTCGGCGCCAGGAGGGTGCAGGA
175▶IleHisAlaSerProAlaGlyThrTyrAlaHisThrValAsnArgAsnTrpTyrSerAspAlaAspValProAlaSerAlaArgGlnGluGlyCysGlnAs
1201 CATCGCTACCGCAGCTCATCTCAACATGGACATTGATGTATCCTGGTGGAGGCGGAAAGTACATGTTTCGCATGGGAACCCAGACCTGAGTACCA
208▶pIleAlaThrGlnLeuIleSerAsnMetAspIleAspValIleLeuGlyGlyGlyArgLysTyrMetPheArgMetGlyThrProAspProGluTyrPro
1301 GATGACTACAGCAAGTGGGACAGGCTGGACGGGAAGATCTGGTGCAGGAATGGCTGGCGAAGCGCCAGGTCGCCGATGTGTGAACCGCACTG
242▶AspAspTyrSerGlnGlyYThrArgLeuAspGlyLysAsnLeuValGlnGluTyrLeuAlaLysArgGlnGlyAlaArgTyrValTyrAsnArgThrG
1401 AGCTCATGACGGCTTCCCTGGACCCGCTGTGACCCATCTCATGGTCTCTTTGAGCCTGGAGACATGAAATACGAGATCCACCGAGACTCCACACTGGA
275▶IleLeuMetGlnAlaSerLeuAspProSerValThrHisLeuMetGlyLeuPheGluProGlyAspMetLysTyrGluIleHisArgAspSerThrLeuAs
1501 CCCCTCCCTGATGGAGATGACAGAGGCTGCCCTGCCCTGCTGAGCAGGAACCCCGCGCTTCTTCTTCTCGTGGAGGGTGGTGCATCGACCCAGCT
308▶pProSerLeuMetGluMetThrGluAlaAlaLeuArgLeuLeuSerArgAsnProArgGlyPhePheLeuPheValGluGlyYArgIleAspHisGly
1601 CATCCAGAAAGCAGGGCTTACCGGGCACTGACTGAGACGATCATGTCGACGACCCATTGAGAGGGCGGGCCAGCTCACCGAGGAGGACACGCTGA
342▶HisHisGlySerArgAlaTyrArgAlaLeuThrGluThrIleMetPheAspAspAlaIleGluArgAlaGlyGlnLeuThrSerGluGluAspThrLeuS
1701 GCCTGCTACCTGCCAGCACTCCACGCTTCTCCTCGGAGGCTACCCCTCGGAGGGAGCTCCATCTTCCGGCTGGCCCTGGCAAGCCCGGACAG
375▶erLeuValThrAlaAspHisSerHisValPheSerPheGlyYThrProLeuArgGlySerSerIlePheGlyLeuAlaProGlyLeuAlaArgAspAr

BamHI (804)
801 CCAAGTCTTAAAGGGCAGAAGAAGGACAACTGGGGCTGAGATACCCCTGGCTATGGACCGCTTCCCATATGTGGCTGTCCAAGACATACAATGT
75▶IaArgIleLeuLysGlyGlnLysLysAspLysLeuGlyProGluIleProLeuAlaMetAspArgPheProTyrValAlaLeuSerLysThrTyrAsnVal
901 AGACAAATATGTCCAGACAGTGGAGCCACAGCCACGCGCTACCTGTGCGGGTCAAGGGCACTTCCAGACCATTTGGCTTGTGAGTGCAGCGCCCGCTTT
108▶IAspLysHisValProAspSerGlyAlaThrAlaThrAlaTyrLeuLysGlyValLysGlyAsnPheGlnThrIleGlyLeuSerAlaAlaAlaArgPhe
1001 AACAGTGAACACGACAGCGCGCAACGAGGTCTCTCGTGTGAATCGGCAAGAAAGCAGGGAAGTCACTGGGAGTGGTAACCACACACGAGTGC
142▶AsnGlnCysAsnThrThrArgGlyAsnGluValIleSerValMetAsnArgAlaLysLysAlaGlyLysSerValGlyValValThrThrArgValG
1101 AGCACGCTCGCCAGCCACCTACGCCACAGGTAACCGCACTGACTCGGACCGCCGCTTTCGAGCTGGAGCCGACGTGCTCGGCGCCAGGAGGGTGCAGGA
175▶IleHisAlaSerProAlaGlyThrTyrAlaHisThrValAsnArgAsnTrpTyrSerAspAlaAspValProAlaSerAlaArgGlnGluGlyCysGlnAs
1201 CATCGCTACCGCAGCTCATCTCAACATGGACATTGATGTATCCTGGTGGAGGCGGAAAGTACATGTTTCGCATGGGAACCCAGACCTGAGTACCA
208▶pIleAlaThrGlnLeuIleSerAsnMetAspIleAspValIleLeuGlyGlyGlyArgLysTyrMetPheArgMetGlyThrProAspProGluTyrPro
1301 GATGACTACAGCAAGTGGGACAGGCTGGACGGGAAGATCTGGTGCAGGAATGGCTGGCGAAGCGCCAGGTCGCCGATGTGTGAACCGCACTG
242▶AspAspTyrSerGlnGlyYThrArgLeuAspGlyLysAsnLeuValGlnGluTyrLeuAlaLysArgGlnGlyAlaArgTyrValTyrAsnArgThrG
1401 AGCTCATGACGGCTTCCCTGGACCCGCTGTGACCCATCTCATGGTCTCTTTGAGCCTGGAGACATGAAATACGAGATCCACCGAGACTCCACACTGGA
275▶IleLeuMetGlnAlaSerLeuAspProSerValThrHisLeuMetGlyLeuPheGluProGlyAspMetLysTyrGluIleHisArgAspSerThrLeuAs
1501 CCCCTCCCTGATGGAGATGACAGAGGCTGCCCTGCCCTGCTGAGCAGGAACCCCGCGCTTCTTCTTCTCGTGGAGGGTGGTGCATCGACCCAGCT
308▶pProSerLeuMetGluMetThrGluAlaAlaLeuArgLeuLeuSerArgAsnProArgGlyPhePheLeuPheValGluGlyYArgIleAspHisGly
1601 CATCCAGAAAGCAGGGCTTACCGGGCACTGACTGAGACGATCATGTCGACGACCCATTGAGAGGGCGGGCCAGCTCACCGAGGAGGACACGCTGA
342▶HisHisGlySerArgAlaTyrArgAlaLeuThrGluThrIleMetPheAspAspAlaIleGluArgAlaGlyGlnLeuThrSerGluGluAspThrLeuS
1701 GCCTGCTACCTGCCAGCACTCCACGCTTCTCCTCGGAGGCTACCCCTCGGAGGGAGCTCCATCTTCCGGCTGGCCCTGGCAAGCCCGGACAG
375▶erLeuValThrAlaAspHisSerHisValPheSerPheGlyYThrProLeuArgGlySerSerIlePheGlyLeuAlaProGlyLeuAlaArgAspAr

StuI (1805)
1801 GAAGGCTACACGGTCTCTCTATACGGAACGGTCCAGGCTATGTGCTCAAGGACGGCGCCCGGGATGTTACCGAGAGCGAGCGGGAGCCCGAG
408▶GlyAlaTyrThrValLeuLeuTyrGlyAsnGlyProGlyTyrValLeuLysAspGlyAlaArgProAspValThrGluSerGluSerGlySerProGlu
1901 TATCGGACAGCTGACAGTGCAGGCTGACGAGAGAGACCCAGCGAGGAGCAGTGGCGGGTGTTCGCGCGCCGCGCCAGCTGGTTCACGGCG
442▶TyrArgGlnGlnSerAlaValProLeuAspGlyGluThrHisAlaGlyGluAspValAlaValPheAlaArgGlyProGlnAlaHisLeuValHisGlyY
2001 TGCAGGACAGACCTTCATAGCGCAGTCAATGGCTTCCGCGCTGCTGGAGCCCTACACCGCTGCGACCTGCGGCCCGCCCGCCAGCACCAGCAGC
475▶AlGlnGluGlnThrPheIleAlaHisValMetAlaPheAlaAlaCysLeuGluProTyrThrAlaCysAspLeuAlaProProAlaGlyThrThrAspAl

Eco147I (1805)
1801 GAAGGCTACACGGTCTCTCTATACGGAACGGTCCAGGCTATGTGCTCAAGGACGGCGCCCGGGATGTTACCGAGAGCGAGCGGGAGCCCGAG
408▶GlyAlaTyrThrValLeuLeuTyrGlyAsnGlyProGlyTyrValLeuLysAspGlyAlaArgProAspValThrGluSerGluSerGlySerProGlu
1901 TATCGGACAGCTGACAGTGCAGGCTGACGAGAGAGACCCAGCGAGGAGCAGTGGCGGGTGTTCGCGCGCCGCGCCAGCTGGTTCACGGCG
442▶TyrArgGlnGlnSerAlaValProLeuAspGlyGluThrHisAlaGlyGluAspValAlaValPheAlaArgGlyProGlnAlaHisLeuValHisGlyY
2001 TGCAGGACAGACCTTCATAGCGCAGTCAATGGCTTCCGCGCTGCTGGAGCCCTACACCGCTGCGACCTGCGGCCCGCCCGCCAGCACCAGCAGC
475▶AlGlnGluGlnThrPheIleAlaHisValMetAlaPheAlaAlaCysLeuGluProTyrThrAlaCysAspLeuAlaProProAlaGlyThrThrAspAl

BglIII (2138) BsrGI (2162)
2101 CGCGCACCCGGCGGCTCCCGTCCCAAGCGTCTGGATAGATCTGGTGTAAAGCCTTGCATATGTACAGTCCCAGAAGTATCATCTGCTTCTATCTCC
508▶AlaHisProGlyYArgSerArgSerLysArgLeuAspArgSerGlyCysLysProCysIleCysThrValProGluValSerSerValPheIlePhePro
1▶GlyCysLysProCysIleCysThrValProGluValSerSerValPheIlePhePro

BbrPI (2247)
2201 CCAAAGCCCAAGGATGTCTCACCATTACTGACTCCTAAGGTACAGTGTGTGGTGGTAGACATCAGCAAGGATGATCCCGAGGTCAGTTCAGTGGT
542▶ProLysProLysAspValLeuThrIleThrLeuThrProLysValThrCysValValAlaAspIleSerLysAspAspProGluValGlnPheSerTrpP
20▶ProLysProLysAspValLeuThrIleThrLeuThrProLysValThrCysValValAlaAspIleSerLysAspAspProGluValGlnPheSerTrpP
2301 TTGTAGATGATGTGGAGTGCACACAGCTCAGACGCAACCCGGGAGGAGCAGTTCACAGCACTTCCGCTCAGTCACTGAACTTCCATCATGCACCA
575▶heValAspAspValGluValHisThrAlaGlnThrGlnProArgGlyGluGlnPheAsnSerThrPheArgSerValSerGluLeuProIleMetHisGly
53▶heValAspAspValGluValHisThrAlaGlnThrGlnProArgGlyGluGlnPheAsnSerThrPheArgSerValSerGluLeuProIleMetHisGly
2401 GGACTGGCTCAATGGCAAGGAGTTCAAATGCAGGTTCAACAGTGCAGCTTCCCTGCCCCATCGAGAAAACCATCTCCAAAACCAAAGGACAGCCGAAG
608▶nAspTrpLeuAsnGlyLysGluPheLysCysArgValAsnSerAlaAlaPheProAlaProIleGluLysThrIleSerLysThrLysGlyYArgProLys
86▶nAspTrpLeuAsnGlyLysGluPheLysCysArgValAsnSerAlaAlaPheProAlaProIleGluLysThrIleSerLysThrLysGlyYArgProLys

BsrGI (2511)
2501 GCTCCACAGGTGACACCATCCACCTCCCAAGGAGCAGATGGCCAAGGATAAAGTCACTGACCTGCATGATAACAGACTTCTTCCCTGAAGACATTA
642▶AlaProGlnValTyrThrIleProProProLysGluGlnMetAlaLysAspLysValSerLeuThrCysMetIleThrAspPhePheProGluAspIleT
120▶AlaProGlnValTyrThrIleProProProLysGluGlnMetAlaLysAspLysValSerLeuThrCysMetIleThrAspPhePheProGluAspIleT
2601 CTGTGGAGTGGCAGTGGAAATGGGACGCCAGCGGAGAAGTACAAGAACACTCAGCCCATCATGGACACAGATGGCTTCTACTTCTGCTACAGCAAGCTCAA
675▶hrValGluTyrGlnTrpAsnGlyGlnProAlaGluAsnTyrLysAsnThrGlnProIleMetAspThrAspGlySerTyrPheValTyrSerLysLeuAs
153▶hrValGluTyrGlnTrpAsnGlyGlnProAlaGluAsnTyrLysAsnThrGlnProIleMetAspThrAspGlySerTyrPheValTyrSerLysLeuAs
2701 TGTGCAGAAGAGCACTGGGAGGACGAAATACTTCCCTGCTGTTGATACATGAGGGCTGCACAACCACTGAGAGGAGCTTCCACTCTCCACTCT
708▶nValGlnLysSerAsnTrpGluAlaGlyAsnThrPheThrCysSerValLeuHisGluGlyLeuHisAsnHisHisThrGluLysSerLeuSerHisSer
186▶nValGlnLysSerAsnTrpGluAlaGlyAsnThrPheThrCysSerValLeuHisGluGlyLeuHisAsnHisHisThrGluLysSerLeuSerHisSer

NheI (2823)
2801 CCTGGTAAATGATCCAGTGCCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTT
742▶ProGlyLys...
220▶ProGlyLys...

HpaI (2963)
2901 ATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTAAGCTGCAATAAACAAGTTAAACAACAACAATTGCATTCATTTTATGTTTCAGTTT

AseI (3060)
3001 AGGGGGAGGTGGGGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTAATTTCTAAATACAGCATAGCAAACTTTAACTCCAAA

3101 TCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTTC

3201 ATGGAGTTTAAAGATATAGTGATTTTTCCCAAGGTTTGAAGTCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCCACATTCCTTTTTTAGTAA

SwaI (3317)

3301 ATATTCAGAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTT

3401 AGTAGTTGGACTTAGGGAAACAAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCTGCCACAAAGT
125↓●●●AspGlnGluGluAlaValPheHis

3501 GCACGAGTTGCCGGCCGGTTCGCGCAGGGCAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTCATGGCCGGCCGGAGGGCTCCCGAAGTTCTGT
116↓sValCysAsnGlyAlaProAspArgLeuAlaPheGluArgGlyTrpProGlnGluGlyIleGluThrMetAlaProGlySerAlaAspArgPheAsnThr

3601 GGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCCGCGCACCCACCCAGGCCAGGGTGTGTCCGGCACCACTGGTCTGGACCGCGCTG
83↓SerValValGluSerTrpGluAlaTyrLeuGluAspLeuGlyArgValTrpValTrpAlaLeuThrAsnAspProValValGluAspGlnValAlaSerI

SgrAI (3730)

3701 ATGAACAGGGTACGTCGTCGCCGACACACCGCGAAGTCTCTCCACGAAGTCCCGGAGAAACCCGAGCCGGTCCGAGAACTCGACCGTCCGG
49↓IlePheLeuThrValAspAspArgValValGlyAlaPheAspAspGluValPheAspArgSerPheGlyLeuArgAspThrTrpPheGluValAlaGlyAla

3801 CGACGTCGCGCGGTGAGCACCGAACGGCACTGGTCAACTTGGCCATGATGGCTCTCctgtcaggagaggaagagaagaaggttagtacaattgCT
16↓aValAspArgAlaThrLeuValProValAlaSerThrLeuLysAlaMet

AseI (3944)

3901 ATAGTGAGTTGATTATACTATGCGATATACATATGCCAATGATTAATTGTCAAACCTAGGGCTGCAgggttcatagtgccacttttctctgactgcccc

HindIII (4070)

4001 tctcctgccaccctttccaggcatagacagtcagtgacttacCAAACCTCACAGGAGGGAGAAGGCAGAAGCTTGAGACAGACCCGCGGGACCGCGAA

StuI (4172)
Eco147I (4172)

4101 CTGCGAGGGGACGTGGCTAGGGCGGCTTCTTTTATGGTGCGCCGCCCTCGGAGGCAGGGCGCTCGGGGAGGCCTAGCGCCAATCTCGGTTGGCAGGAG

BspEI (4228)

4201 CGCGGGCCGAAGCCGTGCCTGACCAATCCGGAGCACATAGGAGTCTCAGCCCCCGCCCCAAAGCAAGGGGAAGTACACGCGCTGTAGCGCCAGCGTGT

SpeI (4335)

4301 TGTGAAATGGGGCTTGGGGGGTTGGGGCCCTGACTAGTCAAACAACCTCCCAATTGACGTCAATGGGGTGGAGACTTGAAATCCCCTGTAGTCAAA

SnaBI (4465)

4401 CGCTATCCACGCCCAATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGG

4501 TCATGTACTGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGC

4601 AGTTTACCCTAAATACTCCACCAATTGACGTCAATGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCAATTATTGACGTCAATGGGCGGGGCTG

PacI (4758)

SdaI (4751)

4701 TTGGGCGGTACGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCG

4801 TAAAAAGGCCCGTGTGCTGGCGTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCAAACCCGACAGGA

4901 CTATAAGATACCAGCGTTTTCCCTGGAAGCTCCCTCGTGCGCTCTCCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGG

5001 GAAGCGTGGCGTTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGTTCGTTCCGCTCAAGCTGGGCTGTGTGCAGAACCCCGTTACGCC

5101 CGACCGCTGCGCTTATCCGGTAACTATCGTCTTGTAGTCCAAACCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGA

5201 GCGAGGTATGTAGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAG

5301 TTACCTTCGAAAAAGAGTTGGTAGCTTTGATCGGCAAAACAACCCCGTGGTAGCGGTGTTTTTTGTTTGAAGCAGCAGATTACGGCGAGAAA

PacI (5498)

5401 AAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGCTGACGCTCAGTGGAAACGAAACTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTA

SwaI (5506) NotI (5514)

5501 ACATTTAAATCAGCGCCGCAATAAAATATCTTTATTTTATTACATCTGTGTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAA

5601 CAAAACGAAACAAAACAACCTAGCAAAATAGGCTGTCCCAGTGAAGTGCAGGTGCCAGAACATTTCTCTATCGAA