

pFUSE-rIgG-Fc2 (IL2ss)

Plasmid designed for the construction of Fc-Fusion proteins

Catalog # pfuse-rfc2

For research use only

Version 20K05-MM

PRODUCT INFORMATION

Content:

- 20 µg of pFUSE-rIgG-Fc2 (IL2ss) plasmid provided as lyophilized DNA
- 1 ml of Zeocin™ (100 mg/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 and lyophilized.

GENERAL PRODUCT USE

pFUSE-Fc is a family of plasmid developed to facilitate the construction of Fc-Fusion proteins by fusing a sequence encoding a given protein to the Fc region of an immunoglobulin.

pFUSE-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, CHO cells, monkey COS cells and human embryonic kidney (HEK)293 cells. These cells are commonly used in protein purification systems.

pFUSE-Fc2 (IL2ss) plasmids allow the secretion of Fc-Fusion proteins. They contain the IL2 signal sequence (IL2ss) for the generation of Fc-Fusion proteins derived from proteins that are not naturally secreted. As Fc-Fusion proteins are secreted, they can be easily detected in the supernatant of pFUSE-Fc-transfected cells by SDS-PAGE. Furthermore, functional domains can be identified by immunoblotting and ligand blotting.

Fc-Fusion proteins can be easily purified by single-step protein A or protein G affinity chromatography.

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.
- **MCS:** The multiple cloning site contains several restriction sites that are compatible with many other enzymes, thus facilitating cloning.
- **IL2 ss:** The IL2 signal sequence contains 20 amino acids 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 antigenic protein.
- **rIgG Fc (rabbit):** The Fc 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 two parts of the Fc-Fusion protein, 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.

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

Zeocin™ usage

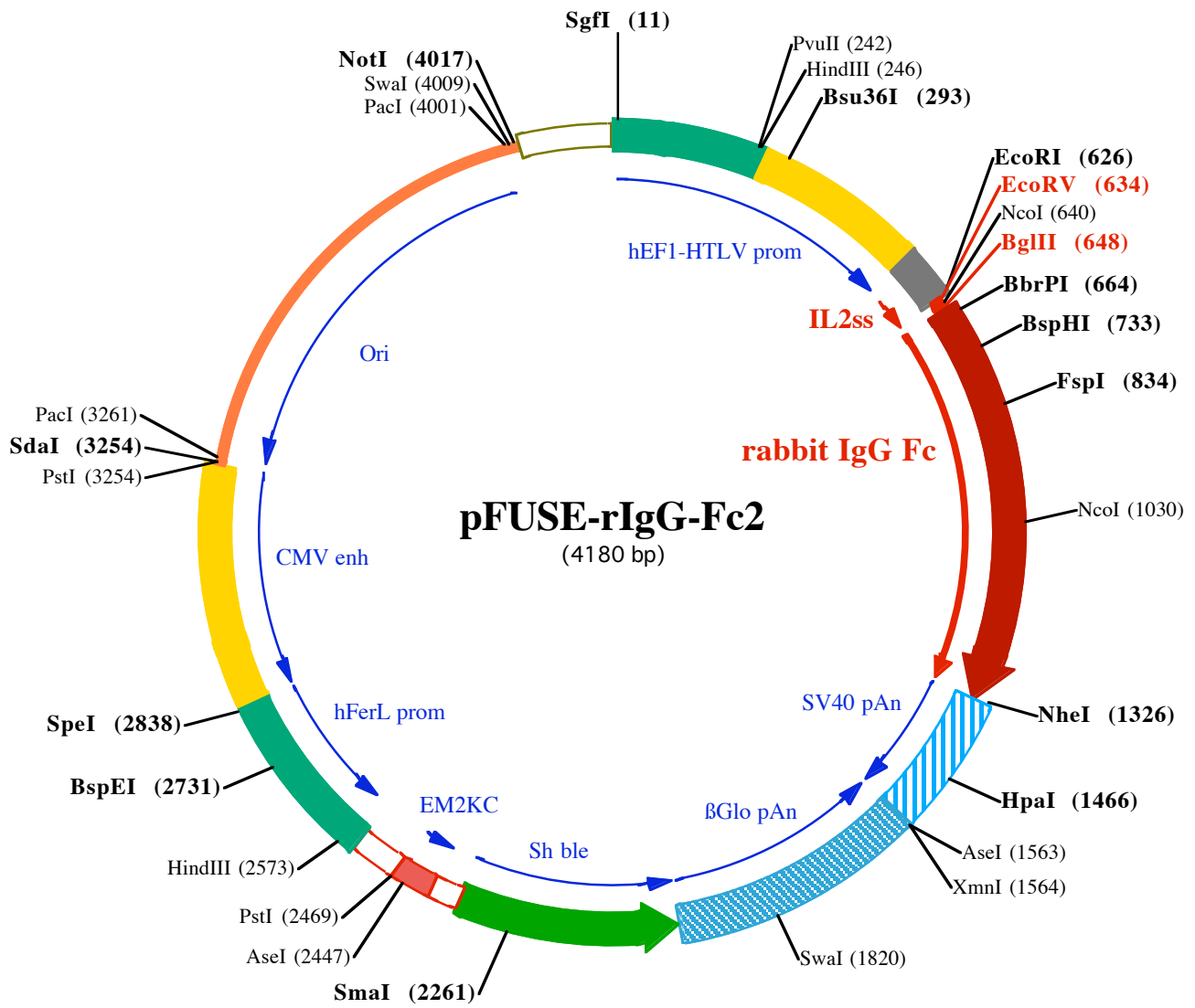
This antibiotic can be used for *E. coli* at 25 µg/ml in liquid or solid media and at 50-200 µg/ml to select Zeocin™-resistant mammalian cells.

References:

1. Kim, D.W. *et al.* (1990). *Gene* 2: 217-223.
2. Takebe, Y. *et al.* (1988). *Mol. Cell Biol.* 1: 466-472.
3. Carswell, S., and Alwine, J.C. (1989). *Mol. Cell Biol.* 10: 4248-4258.
4. Yu J & Russell JE. (2001). *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 Hong Kong: +852 3622-3480
E-mail: info@invivogen.com



SgfI (11)
1 GGATCTGCGATCGCTCCGGTCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGTTCGCAATTGAACGGGTGCCTA
101 GAGAAGGTGGCGGGGTAACCTGGGAAAGTGATGTCGTGTACTGGCTCCGCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (246)
PvuII (242)
Bsu36I (293)
201 GTGAACGTTCTTTTTCGCAACGGGTTTGGCCGAGAAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCCTTACGCGCCCGCCCTACCTGAGGCC
301 GCCATCCACGCGGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCTGAATCGCTCCGCGCTTAGGTAAGTTTAAAGCTCAGGTCGAGACC
401 GGGCCTTTGTCGGCGCTCCCTTGAGCCTACCTAGACTCAGCGGGCTCCACGCTTTGCTGACCTGCTTGTCTCAACTCTACGCTTTGTTTCGTTT
501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGGCCCTACCTGAGATCAcggcGAAGGAGGGCCACCATGTACAGGATGCAACTCCTGTCTTGCA
1►MetTyrArgMetGlnLeuLeuSer CysI

EcoRV (634)
EcoRI (626) **NcoI (640)** **BglIII (648)** **BbrPI (664)**
601 TTGCACTAAGTCTTGCACCTGTACGAATTCGATATCGGCCATGGTTAGATCTAGCAAGCCACGTGCCACCCCTGAATCTTGGGGGACCGTCTGT
10►IleAlaLeuSerLeuAlaLeuValThrAsnSer 1►SerLysProThrCysProProRoGluLeuLeuGlyProSerVa

BspHI (733)
701 CTTTCATCTTCCCAAAACCAAGGACACCTCATGATCTCAGCACCCCGAGGTCACATGCGTGGTGGACGTGAGCGAGGATGACCCCGAGGTC
16►IPheIlePheProProLysProLysAspThrLeuMetIleSerArgThrProGluValThrCysValValValAspValSerGlnAspAspProGluVal

FspI (834)
801 CAGTTCACATGGTACATAAACAACGAGCAGTGGCGCACCCTCCCGCGCGCTACGGGAGCAGAGTTCAACAGCAGATCCGCGTGGTCAGCACCTCC
50►GlnPheThrTrpTyrIleAsnAsnGluGlnValArgThrAlaArgProProLeuArgGluGlnGlnPheAsnSerThrIleArgValValSerThrLeuP
901 CCATCGCGCACAGGACTGGCTGAGGGCAAGGAGTTCAAGTGCAAAGTCCACAACAAGGCACTCCCGCCCCATCGAGAAAACCATCTCCAAAGCCAG
83►roIleAlaHisGlnAspTrpLeuArgGlyLysGluPheLysCysLysValHisAsnLysAlaLeuProAlaProIleGluLysThrIleSerLysAlaAr

NcoI (1030)
1001 AGGGCAGCCCTGGAGCCGAAGGTCTACACCATGGGCCCTCCCGGGAGAGTGGAGCAGCAGGTCGGTCACTGACCTGCATGATCAACGGCTTCTAC
116►gGlyGlnProLeuGluProLysValTyrThrMetGlyProProArgGluGluLeuSerSerArgSerValSerLeuThrCysMetIleAsnGlyPheTyr
1101 CCTTCCGACATCTCGTGGAGTGGGAGAAGAAGCGGAGGCAAGTCAAGACACCGCGCGCTGCTGGACAGCGACGGCTCCTACTTCTCT
150►ProSerAspIleSerValGluTrpGluLysAsnGlyLysAlaGluAspAsnTyrLysThrThrProAlaValLeuAspSerAspGlySerTyrPheLeuT
1201 ACAGCAAGCTCTCAGTCCACGAGTGGTGGCAGCGGGGCGAGCTTCCACCTGCTCCGTGATGCACGAGGCTTGCACAACCACTACACGCAAGATC
183►yrserLysLeuSerValProThrSerGluTrpGluArgGlyAspValPheThrCysSerValMetHisGluAlaLeuHisAsnHisTyrThrGlnLysSe

NheI (1326)
1301 CATCTCCGCTCTCCGGTAAATGAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAAGTGAATGCAATGAAAAAATGC
216►rIleSerArgSerProGlyLys•••

HpaI (1466)
1401 TTTATTTGTAAATTTGTGATGCTATTGCTTTATTTGTAACCTATAAGCTGCAATAAACAAGTTAAACAACAATTCATTCTTTATGTTTCAGG

AseI (1563)
XmnI (1564)
1501 TTCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATTAATCTAAAATACAGCATAGCAAACTTTAACCTCC
1601 AAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTCT
1701 TTCATGGAGTTAAGATATAGTGATTTTCCCAAGGTTTGAAGTACTCTTCAATTTCTTATGTTTTAAATGCAGTACCTCCACATCCCTTTTATG

SwaI (1820)
1801 AAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAG
1901 TTTAGTAGTTGGACTTAGGAAACAAAGGAACCTTTAATAGAATTTGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTGCCACAA
125►•••AspGlnGluGluAlaValPh

SmaI (2261)
2201 AGTGCACGAGTTGCGGCCGGTTCGCGCAGGCGCAACTCCCGCCCAACGCTGCTCGCGATCTCGGTATGGCCGGCGGGGCTCCCGAAGTT
117►eHisValCysAsnGlyAlaProAspArgLeuAlaPheGluArgGlyTrpProGlnGlyIleGluThrMetAlaProGlySerAlaAspArgPheAsn
2101 CGTGGACACGACTCGGACACTCGGCGTACAGCTCGTCCAGGCGCGCACCCACCCAGGCCAGGGTGTGTCGGCACCACTGCTGCTGGACCGCG
84►ThrSerValValGluSerTrpGluAlaTyrLeuGluAspLeuGlyArgValTrpValIleuThrAsnAspProValValGlnAspGlnValAlaIAs

AseI (2447) **PstI (2469)**
2401 gCTATAGTGAGTTGATTATACTATGCAGATATACTATGCCAATGATTAATTGTCAAACCTAGGGCTGCAgggttcatagtgccacttttctcgtcactgcc
17►yAlaValAspArgAlaThrLeuValProValAlaSerThrLeuLysAlaMet

HindIII (2573)
2501 ccatctctctgccaccctttccaggcatagacagtcagtgacttacCAAACCTACAGGAGGAGAAGGCAGAAGCTTGAGACAGACCCCGGGACCGCC
2601 GAACTGCGAGGGGACGTGGCTAGGGGGCTTTTATGTTGCGCGGCCCTCGGAGGCAAGGGCTCGGGGAGGCTAGCGCCCAATCTGCGGTGGCAG

BspEI (2731)
2701 GAGGCGGGCCGAAGCCGTGCCTGACCAATCCGGAGCACATAGGAGTCTCAGCCCCCGCCCAAGCAAGGGGAAGTACGCGCTGTAGCGCCAGCG

SpeI (2838)
2801 TGTTGTAAATGGGGCTTGGGGGGTGGGGCCCTGACTAGTCAAACAAACTCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCGTGAGTCA
2901 AACCGCTATCCACGCCATTGATGTAAGTCCAAAACCGCATCATGGTAAATAGCGATGACTAATACGTAGATGTAAGTCCAAAGTGGAAAGTCCATA
3001 AGGTCATGTAAGTGGCATAATGCCAGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGTACTGCAAGTGG
3101 GGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGAAACATACGTCAATTTGACGTCAATGGCGGGGG

PacI (3261)

PstI (3254)

SdaI (3254)

3201 TCGTTGGGCGGT CAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAA
3301 CCGTAAAAAGGCCGCTTGTGGCGTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACA AAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACA
3401 GGA CTATAAAGATACCAGGCGTTTTCCCTGGAAGCTCCCTGTGCGCTCTCCTGTTCCGACCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTT
3501 CGGGAAGCGTGGCGCTTTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCA
3601 GCCCGACCGCTGCGCTTATCCGTA ACTATCGTCTTGAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGC
3701 AGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGC
3801 CAGTTACCTTCGGAAAAGAGTTGGTAGCTTTGATCCGGCAAACAACCCCGCTGGTAGCGGTGGTTTTTTTGTTCGAAGCAGCAGATTACGCGCAG

PacI (4001)

3901 AAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAA

SwaI (4009)

NotI (4017)

4001 TTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTAATCGTAACTAACATACGCTCTCCATCA
4101 AAACAAAACGAAACAAAACAAACTAGCAAATAGGCTGTCCCGAGTGAAGTGCAGGTGCCAGAACATTTCTCTATCGAA