

# pFUSE-mIgG3-Fc2

Plasmid containing a mouse IgG3 Fc region

Catalog # pfuse-mg3fc2

For research use only

Version 20K05-MM

## PRODUCT INFORMATION

### Content:

- 20 µg of pFUSE-mIgG3-Fc2 (ssIL2) 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 the effector region of a protein to the Fc region of an immunoglobulin G (IgG).

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, cells that 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.

InvivoGen provides pFUSE-Fc vectors featuring Fc regions from different species and isotypes. Three murine isotypes are available: IgG1, IgG2a and IgG3. The Fc region mediates effector functions, such as antibody-dependent cellular cytotoxicity (ADCC) and complement-dependent cytotoxicity (CDC). In ADCC, the Fc region of an antibody binds to Fc receptors (FcγRs) on the surface of immune effector cells such as natural killers and macrophages, leading to the phagocytosis or lysis of the targeted cells. In CDC, the antibodies kill the targeted cells by triggering the complement cascade at the cell surface. IgG isoforms exert different levels of effector functions increasing in the order of mIgG1<mIgG3<mIgG2a.

## PLASMID FEATURES

- **mIgG3 Fc (mouse):** 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. The Fc region of mouse IgG3 mediates high ADCC and low CDC<sup>1</sup>.
- **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1α (EF-1α) core promoter<sup>2</sup> and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat<sup>3</sup>. 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.
- **IL2 ss:** The IL2 signal sequence contains 22 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.
- **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 mRNA<sup>4</sup>.
- **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 transcription<sup>5</sup>.

### References:

1. Dangel JL. *et al.*, Segmental flexibility and complement fixation of genetically engineered chimeric human, rabbit and mouse antibodies. *EMBO J.* 7(7):1989-94.
2. Kim DW *et al.* 1990. Use of the human elongation factor 1 alpha promoter as a versatile and efficient expression system. *91(2):217-23.*
3. 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.
4. 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.
5. 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 Hong Kong: +852 3622-3480  
E-mail: [info@invivogen.com](mailto: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 H<sub>2</sub>O. 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.

## RELATED PRODUCTS

Product	Catalog Code
Zeocin™	ant-zn-1

---

### TECHNICAL SUPPORT

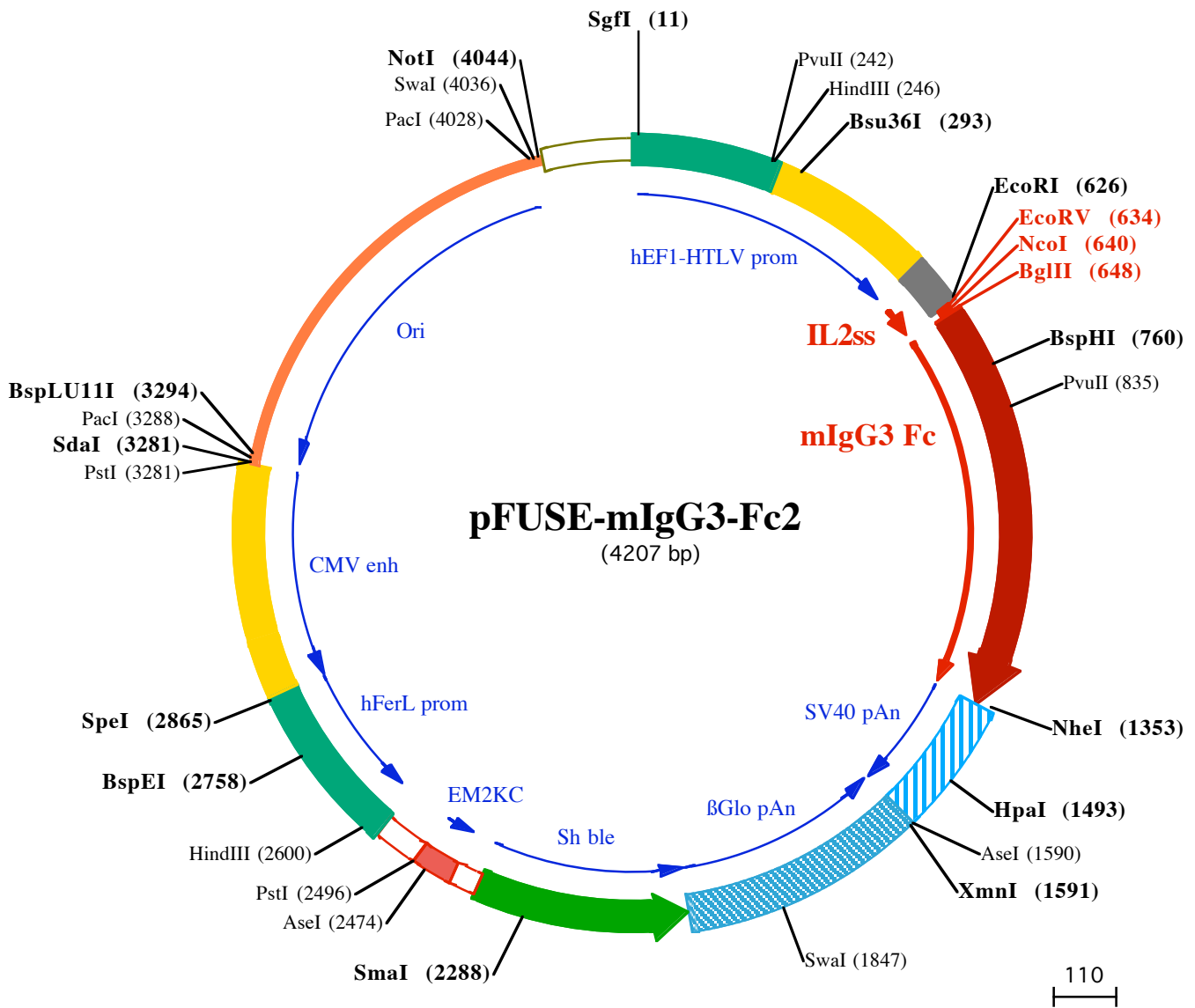
InvivoGen USA (Toll-Free): 888-457-5873

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](mailto:info@invivogen.com)



**SgfI (11)**  
1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGCAGAGCGCACATCGCCACAGTCCCGAGAAGTTGGGGGAGGGGTGCGCAATTGAACGGGTGCCTA  
101 GAGAAGGTGGCGCGGGGTAACCTGGGAAAGTGATGCTGTACTGGCTCCGCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

**HindIII (246)**  
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACAACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTTCACGCGCCCGCCCTACCTGAGGCC  
**PvuII (242)** **Bsu36I (293)**

301 GCCATCCACGCGGTTGAGTCGCGTCTGCCGCTCCCGCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTAAAGCTCAGGTCGAGACC  
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCGGGCTCTCCACGCTTTGCTGACCTGCTTCACTACGTCTTTGTTTCGTTT  
501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTACCAGGCTACCTGAGATCAcggcGAAGGAGGGCCACCATGTACAGGATGCAACTCTGTCTTGCA  
1►MetTyrArgMetGlnLeuLeuSer CysI

**EcoRV (634)** **BglII (648)**  
601 TTGCACTAAGTCTTGCACCTGTGCACGAATTCGATATCGGCCATGGTTAGATCTCTCTAGAATACCCAGCCAGTACCCCGGTTCTTCATGCCACC  
10►LeuAlaLeuSerLeuAlaLeuValThrAsnSer 1►ProArgIleProLysProSerThrProProGlySerSerCysProPro

**EcoRI (626)** **NeoI (640)**  
701 TGGTAACATCTTGGGTGGACCATCCGTCTTCTTCCCCCAAAGCCAAAGGATGCACTCATGATCTCCCTAACCCCAAGGTTACGTGTGGTGGTG  
16►oGlyAsnIleLeuGlyGlyProSerValPheIlePheProProLysProLysAspAlaLeuMetIleSerLeuThrProLysValThrCysValValVal

**BspHI (760)**  
801 GATGTGAGCGAGGATGACCCAGATGTCCATGTGACGTGGTTGTGGACAACAAGAAGTACACACAGCTGGACACAGCCCGTGAAGCTCAGTACAACA  
50►AspValSerGluAspAspProAspValHisValSerTrpPheValAspAsnLysGluValHisThrAlaTrpThrGlnProArgGluAlaGlnTyrAsnS  
901 GTACCTCCGAGTGGTCACTGCCCTCCCATCCAGCACCAGGACTGGATGAGGGCAAGGAGTTCAAATGCAAGGTCAACAACAAGCCCTCCAGCCCC  
83►erThrPheArgValValSerAlaLeuProIleGlnHisGlnAspTrpMetArgGlyLysGluPheLysCysLysValAsnAsnLysAlaLeuProAlaPr  
1001 CATCGAGAGAACCATCTCAAAACCCAAAGGAAGAGCCAGACACCTCAAGTATACACCATACCCCACTCTGTAACAAATGCAAGAAAGGTTAGT  
116►oIleGluuArgThrIleSerLysProLysGlyArgAlaGlnThrProGlnValTyrThrIleProProProArgGluGlnMetSerLysLysLysValSer  
1101 CTGACCTGCCGTGGTCAACCACTTCTTCTGAAGCCATCAGTGTGGAGTGGGAAAGGAACCGGAGAAGTGGAGCAGGATTACAAGAACACTCCACCCATCC  
150►LeuThrCysLeuValThrAsnPhePheSerGluAlaIleSerValGluTrpGluuArgAsnGlyGluLeuGluGlnAspTyrLysAsnThrProProIleLe  
1201 TGGACTCAGATGGGACTACTTCTCTACAGCAAGCTCACTGTGGATACAGACAGTGGTTGCAAGGAGAAATTTTACCTGCTCCGTGGTGCATGAGGC  
183►LeuAspSerAspGlyThrTyrPheLeuTyrSerLysLeuThrValAspThrAspSerTrpLeuGlnGlyIlePheThrCysSerValValHisGluAl

**NheI (1353)**  
1301 TCTCCATAACCACACACAGAAGAACCTGTCTCGCTCCCTGGTAAATGAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAAC  
216►aLeuHisAsnHisHisThrGlnLysAsnLeuSerArgSerProGlyLys•••

**HpaI (1493)**  
1401 ACAACTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGTGAATAAACAAGTTAAACAACA

**AseI (1590)**  
1501 ACAATTGCATTCAATTTATGTTTCAGGTTTCAGGGGAGGTGTGGAGGTTTTTTTAAAGCAAGTAAAACCTCTACAATGTGGTATGGAATTAATCTAAA  
**XmnI (1591)**

1601 ATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTG  
1701 CATTAGCTGTTTGCAGCCTCACCTCTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGGTTTGAAGTCTTTCATTCTTTATGTTTAAATGC

**Swal (1847)**  
1801 ACTGACCTCCACATTCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTTATTAGGCAGAATCCAGATG  
1901 CTCAAGGCCCTTCAATAATCCCCAGTTTAGTAGTTGGACTTAGGAAACAAAGGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTT

2001 ATCCTCAGTCTGCTCTGTCACAAAGTGCACGAGTTGCCGGCGGGTGCAGCAGGGCAACTCCCGCCCAACGGCTGCTCGCCGATCTCGGTGAT  
125►•••AspGlnGluGluAlaValPheHisValCysAsnGlyAlaProAspArgLeuAlaPheGluuArgGlyTrpProGlnGluGlyIleGluuThrMet  
2101 GCGCGCCCGGAGGCTCCCGAAGTTCGTGGACAGCCTCCGACCCTCGGCTGACGCTGCTCAGGCCGCGCACCCACACCAGGCCAGGTTGTTG  
93►AlaProGlySerAlaAspArgPheAsnThrSerValValGluuSerTrpGluAlaTyrLeuGluuAspLeuGlyArgValTrpValTrpAlaLeuThrAsnA

**SmaI (2288)**  
2201 TCCGGCACCACTGGTCTGGACCGCTGATGAACAGGGTACGTCGTCGCGGACCAACCGGCAAGTCTGCTCCACGAAGTCCCGGAGAACCCTCGA  
59►spProValValGluAspGlnValAlaSerIlePheLeuThrValAspAspArgValValGlyAlaPheAspAspGluValPheAspArgSerPheGlyLe  
2301 GCCGTCGTTCCAGAACTCGACCGCTCCGGCAGCTCGCGCGGGTGAACCGGAAACCGCACTGGTCAACTTGGCCATGATGGCTCCTCctgtcaggag  
26►uArgAspThrTrpPheGluValAlaGlyAlaValAspArgAlaThrLeuValProValAlaSerThrLeuLysAlaMet

**AseI (2474)** **PstI (2496)**  
2401 aggaagagaagaaggttagtacaattgCTATAGTGAGTTGATTATACTATGCAGATATACTATGCCAATGATTAATGTCAAACCTAGGGCTCGAaggt

**HindIII (2600)**  
2501 tcatagtccacttttctgactgccccatctctgccccctttccaggcatagacagtcagtgacttacAAACTCAGGAGGGAGAAGGCAGA  
2601 AGCTTGAGACAGACCCCGGGACCGCGAACTGCGAGGGAGCTGGCTAGGGCGGCTTCTTTATGTTGCGCCGCGCTCGGAGGCAGGGCTCGGGGA

**BspEI (2758)**  
2701 GGCCTAGCGGCAATCTGCGGTGGCAGGAGCGGGGCCGAAGCGCTGCCTGACCAATCCGGAGCACATAGGAGTCTCAGCCCCCGCCCAAGCAAGG

**SpeI (2865)**  
2801 GGAAGTACGCGCTGTAGCGCCAGCGTGTGTGAAATGGGGCTTGGGGGTTGGGGCCCTGACTAGTCAAAACAACCTCCATTGACGTCAATGGGG  
2901 TGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCACGCCATTGATGACTGCCAAAACCGCATCATGGTAATAGCGATGACTAATACGTAGA  
3001 TGTACTGCAAGTAGGAAAGTCCATAAGGTCATGACTGGGCATAATGCCAGCGGGCCATTTACCGTCAATTGACGTCAATAGGGGCGTACTTGGCAT  
3101 ATGATACACTTGATGACTGCCAAGTGGGCGAGTTTACCATAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGAAACATA

Pacl (3288)  
PstI (3281) **SdaI (3281)** **BspLU111 (3294)**

3201 GTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTCAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCCTGCAGGTTAATTAAGAACATGTG  
3301 AGCAAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGC  
3401 TCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCCTGCCGTTA  
3501 CCGGATACCTGTCCGCCTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCT  
3601 GGGCTGTGTGCACGAACCCCGTTAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTG  
3701 GCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAG  
3801 TATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTTTGATCCGGCAAACAAACCACCGTGGTAGCGGTGGTTTTTT  
3901 TGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTACGGGGTCTGACGCTCAGTGAACGAAAACTCAGT  
Pacl (4028) SwaI (4036) **NotI (4044)**  
4001 TAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGA  
4101 ATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAAATAGGCTGTCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTC  
4201 TATCGAA