

pFUSE-mIgG1-Fc2

Plasmid containing a mouse IgG2 Fc region and the IL-2 signal sequence

Catalog code: pfuse-mg1fc2

<https://www.invivogen.com/pfuse-migg1-fc>

For research use only

Version 20K05-MM

PRODUCT INFORMATION

Contents:

- 20 µg of pFUSE-mIgG1-Fc2 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.

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 that are commonly used in protein purification systems.

pFUSE-Fc plasmids allow the secretion of Fc-Fusion proteins. 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.

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

PLASMID FEATURES

- **mIgG1 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 IgG1 mediates low CDC and no ADCC¹.
- **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1α (EF-1α) core promoter² and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat³. 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.
- **IL-2 ss:** The IL2 signal sequence contains 21 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 mRNA⁴.
- **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 transcription⁵.

1. Kipps TJ. *et al.*, 1985. Importance of immunoglobulin isotype in human antibody-dependent, cell-mediated cytotoxicity directed by murine monoclonal antibodies. *J Exp Med.* 161(1):1-17. 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



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.

RELATED PRODUCTS

Product	Catalog Code
Zeocin™	ant-zn-1

TECHNICAL SUPPORT

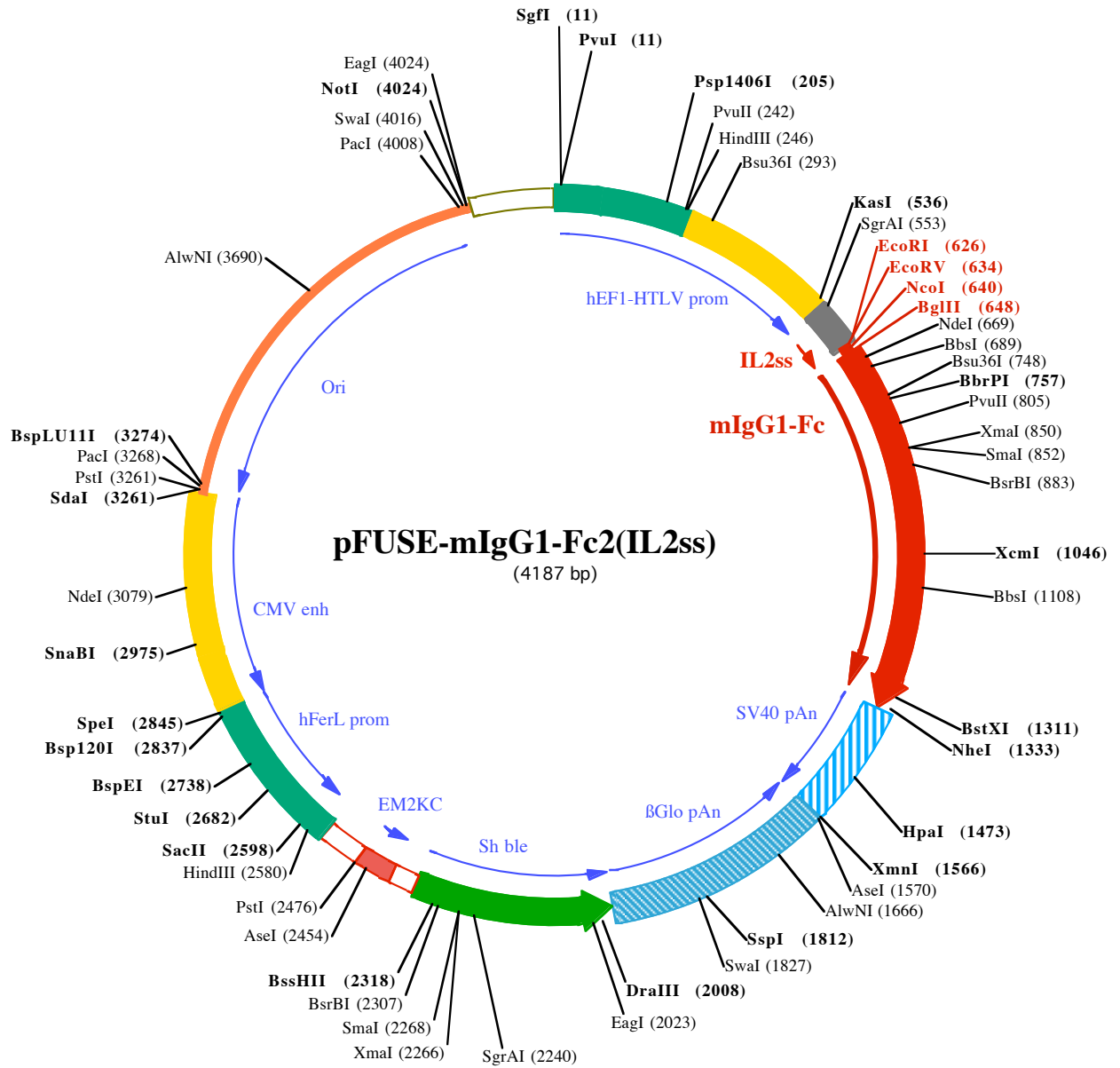
InvivoGen USA (Toll-Free): 888-457-5873

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



100

PvuI (11)
SgfI (11)
 1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
 101 GAGAAAGTGGCGGGGTAAGTGGAAAGTGATGTCGTGACTGGCTCCGCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (246)
Psp1406I (205) **PvuII (242)** **Bsu36I (293)**
 201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACCGCGCCGCCGCCCTACCTGAGGGCC
 301 GCCATCCACGCGGTTGAGTGCAGTCTGCCGCTCCCGCTGTGGTGCTCCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC
 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTCTCAACTCTACGTCTTTGTTTCGTTT

KasI (536) **SgrAI (553)**
 501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCGCTACTCTGAGATCAccggcGAAGGAGGGCCACCATGTACAGGATGCAACTCCTGTCTTGCA
 1►MetTyrArgMetGlnLeuLeuSer CysI

EcoRV (634) **BglIII (648)**
EcoRI (626) **NcoI (640)** **NdeI (669)** **BbsI (689)**
 601 TTGCACTAAGTCTTGCACTTGTCACGAATTCGATATCGGCCATGGTTAGACTCTGGTTGAAGCCTTGCATATGTACAGTCCCAGAAAGTATCATCTGTCTT
 1►IeAlaLeuSerLeuAlaLeuValThrAsnSer 1►GlyCysLysProCysIleCysThrValIProGluValSerSerValPh

BbrPI (757)
Bsu36I (748)
 701 CATCTCCCCCAAAGCCCAAGGATGTGCTCACCATTCTGACTCCTAAGGTCACGTGTGTTGTGGTAGACATCAGCAAGGATGATCCCGAGGTCCAG
 1►eIlePheProProLysProLysAspValLeuThrIleThrLeuThrProLysValThrCysValValValAspIleSerLysAspAspProGluValGln

XmaI (850) **SmaI (852)** **BsrBI (883)**
 801 TTCAGCTGGTTGTAGATGATGTGGAGGTGCACACAGCTCAGACGCAACCCCGGAGGAGCAGTTCAACAGCACTTCCGCTCAGTCAAGTCCCA
 50►PheSerTrpPheValAspAspValGluValHisThrAlaGlnThrGlnProArgGluGluGlnPheAsnSerThrPheArgSerValSerGluLeuProI
 901 TCATGCACCAGACTGGCTCAATGGCAAGGAGTCAAATGCAGGGTCAACAGTGCAGCTTCCCTGCCCATCGAGAAAACCATCTCCAAAACCAAAGG
 83►IleMethisGlnAspTrpLeuAsnGlyLysGluPheLysCysArgValAsnSerAlaAlaPheProAlaProIleGluLysThrIleSerLysThrLysGln

XcmI (1046)
 1001 CAGACCGAAGGCTCCACAGGTGTACACCATTCCACCTCCCAAGGAGCAGATGGCCAAGGATAAAGTCAGTCTGACCTGCATGATAACAGACTTCTCCCT
 116►YargProLysAlaProGluValTyrThrIleProProProLysGluGlnMetAlaLysAspLysValSerLeuThrCysMetIleThrAspPhePhePro

BbsI (1108)
 1101 GAAGACATTACTGTGGAGTGGCAGTGGAAATGGGACGCCAGCGGAGAACTACAAGAACTCAGCCATCATGGACACAGATGGCTCTTACTTCGTCTACA
 150►GluAspIleThrValGluTrpGlnTrpAsnGlyGlnProAlaGluAsnTyrLysAsnThrGlnProIleMetAspThrAspGlySerTyrPheValTyrS

1201 GCAAGCTCAATGTGAGAAGAGCAACTGGGAGCGAGAAATACTTTCACCTGCTCTGTGTACATGAGGGCTGCACAACCACCACTAGAGAAGAGCCT
 183►erLysLeuAsnValGlnLysSerAsnTrpGluAlaGlyAsnThrPheThrCysSerValLeuHisGluGlyLeuHisAsnHisHisThrGluLysSerLe

BstXI (1311) **NheI (1333)**
 1301 CTCCCCTCTCTGGTAAATGATCCAGTGTGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGACAAACCACAACCTAGAATGCAGTGAAA
 216►uSerHisSerProGlyLys•••

HpaI (1473)
 1401 AAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACCAAGTTAAACAACAATTGCATTATTATG

AseI (1570) **XmnI (1566)**
 1501 TTTCAGTTTCAGGGGAGGTGTGGAGGTTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTAATCTAAAATACAGCATAGCAAACTTT

AlwNI (1666)
 1601 AACCTCCAAATCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTTTGCCAATGTGCATTAGCTGTTTGCAGCCTC
 1701 ACCTTCTTTTATGGAGTTAAGATATAGTGTATTTTCCCAAGGTTTGAAGTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCT

SspI (1812) **SwaI (1827)**
 1801 TTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATAT
 1901 CCCCAGTTTAGTGTGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTCT
 125►•••AspGlnGluGlu

DraIII (2008) **EagI (2023)**
 2001 GCCCAAAGTGCACGAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCCCGCCCCAGGCTGCTCGCCGATCTCGGTCATGGCCGGCCCGAGGCGTCCC
 119►IaValPheHisValCysAsnGlyAlaProAspArgLeuAlaPheGluArgGlyTrpProGlnGluGlyIleGluuThrMetAlaProGlySerAlaAspAr

2101 GGAAGTTCGTGGACACGACTCCGACCCTCGCGTACAGCTCGTCCAGGCGCGCACCCACACCCAGGCCAGGGTGTGTCCGGCACCACTGGTCTG
 86►gPheAsnThrSerValValGluSerTrpGluAlaTyrLeuGluAspLeuGlyArgValTrpValTrpAlaLeuThrAsnAspProValValGlnAspGln

SgrAI (2240) **XmaI (2266)** **SmaI (2268)**
 2201 GACCGCTGATGAACAGGGTCACGTCGTCGCCGACCAACCGGCAAGTCTCTCCACGAAGTCCCGGGAGAACCCGAGCCGGTCCGAGAACTCG
 53►ValAlaSerIlePheLeuThrValAspAspArgValValGlyAlaPheAspAspGluValPheAspArgSerPheGlyLeuArgAspThrTrpPheGluuV

BsrBI (2307) **BssHIII (2318)**
 2301 ACCGCTCCGGCGAGCTCGCGCGGTGAGCACCGGAACGGCACTGGTCAACTTGGCCATGATGGCTCCTCctgtcaggagaggaaagagaagaaggttag
 19►aAlaGlyAlaValAspArgAlaThrLeuValProValAlaSerThrLeuLysAlaMet

AseI (2454) **PstI (2476)**
 2401 tacaattgCTATAGTGAGTTGTATTATACTATGCAGATATACTATGCCAATGATTAATTGTCAAACCTAGGGCTGCAGggttcatagtgcacttttctgt

2501 cactgccccatctcctgcccaccctttccaggcatagacagtcagtgacttacCAAACCTCACAGGAGGGAGAAGGCAGAAGCTTGAGACAGACCCCGCGG HindIII (2580) SacII (2598)

2601 GACCGCCGAAGTGCAGGGGACGTGGCTAGGGCGGCTTCTTTTATGGTGCGCCGCCCTCGGAGGCAGGGCGCTCGGGGAGGCCTAGCGGCAATCTGCG StuI (2682)

2701 GTGGCAGGAGCGGGGCCAAGGCCGTGCCTGACCAATCCGGAGCACATAGGAGTCTCAGCCCCCGCCCAAAGCAAGGGGAAGTCACGCGCCTGTAGC BspEI (2738)

2801 GCCAGCGTGTGTGAAATGGGGGCTTGGGGGGTGGGGCCCTGACTAGTCAAAACAAACTCCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCGG SpeI (2845) Bsp120I (2837)

2901 TGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAG SnaBI (2975)

3001 TCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTACCCTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGTGTACTG NdeI (3079)

3101 CCAAGTGGGCAAGTTACCCTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGG

3201 GCGGGGGTCGTTGGGCGGTGAGCCAGGCGGGCCATTTACCCTAAGTTATGTAACGCCTGCAGGTTAATTAAGAACATGTGAGCAAAGCCAGCAAAGG PacI (3268) PstI (3261) SdaI (3261) BspLU11I (3274)

3301 CCAGGAACCGTAAAAAGCCGCTTGTGGCGTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAA

3401 CCCGACAGGACTATAAAGATACCAGCGTTTTCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCGACCCTGCCGCTTACCGGATACCTGTCCGCTTT

3501 CTCCCTTCGGAAGCGTGGCGCTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCC

3601 CCGTTCAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAG AlwNI (3690)

3701 GATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGATTTGGTATCTGCGCTCTG

3801 CTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAACCACCGCTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTA

3901 CGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTACGGGGTCTGACGCTCAGTGGAACGAAAACTCACGTTAAGGGATTTTGGTCATGGC

4001 TAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTTCATTACATCTGTGTGGTTTTTTTTGTGTGAATCGTAACTAACATACGCTC EagI (4024) PacI (4008) SmaI (4016) NotI (4024)

4101 TCCATCAAACAAAACGAAACAAAACAACTAGCAAAATAGGCTGTCCCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA