

pFUSE-mIgG2A-Fc2

Plasmid containing a mouse IgG2A Fc region

Catalog # pfuse-mg2afc2

For research use only

Version 20K05-MM

PRODUCT INFORMATION

Content:

- 20 µg of pFUSE-mIgG2A-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 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

- **mIgG2A 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 IgG2A mediates high ADCC and CDC.
- **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.
- **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.
- **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.
- **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 *Streptallosteichus hindustanus*. The same resistance gene confers selection in both mammalian cells and *E. coli*.
- **BGlo pAn:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription⁴.

References:

1. 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.
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 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

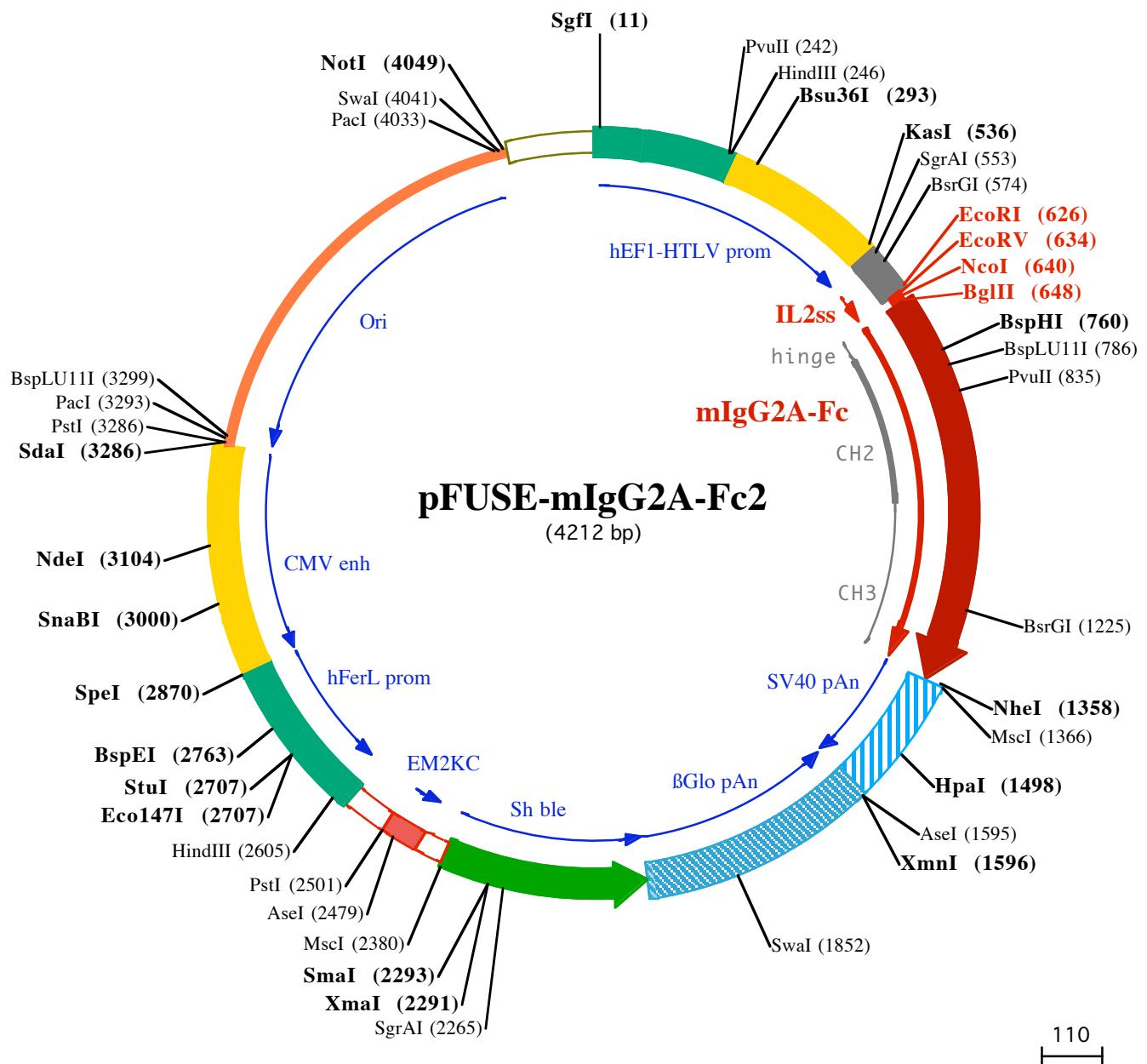
InvivoGen USA (Toll-Free): 888-457-5873

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



SgfI (11)

1 GGATCTGCATCGCCGGTCCCCGTCAAGTGGCAGAGCGCACATGCCACAGTCCCAGAAGTTGGGGGAGGGTCGGCAATTGAACGGTGCCTA

101 GAGAACGGTGGCGCGGGTAAACTGGAAAGTGATCTGTAUTGGCTCGCTTCCGAGGGTGGGGAGAACCGTATAAGTCAGTAGTCGCC

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

201 GTGAACTTCTTTGCACGGGTTGCCAGAACACAGCTAAGCTCGAGGGCTGCATCTCCTCACGGCCCCGCCCTACCTGAGGCC

301 GCCATCCACGCCGGTGAAGCTCGCTGCCCTCCGCTGTGGCTCCTGAACTGCCTCCGTAGGTAAGTTAACGTCAGGTCAGACC

401 GGGCCTTGTCCGGCGTCCCTGGAGCCTACCTAGACTCAGCCGCTCTCACGCTTGCTGACCCGTCACTCTAGTCAGTCTTGTCTT

KasI (536) SgrAI (553) BsrGI (574)

501 TCTGTTCTGCCTTACAGATCCAAGCTGACCGCGCTACCTGAGATCACGGCGAAGGAGGGCCACATGACAGGATGCAACTCCTGCTTGC

1►MetTyrArgMetGl nLeuLeuSer CysI

EcoRV (634) BglII (648)
EcoRI (626) NcoI (640)

601 TTGCACTAAGTCTTGCACTTGCAATTGATATGCCCATGGTAGATCTCCAGAGGGCCACAATCAAGCCCTGCTCTCATGCAAATGCCAGC

10►IeAl aLeuSer LeuAl aLeuVal Thr AsnSer

BspHI (760) BspLU11I (786)

701 ACCTAACCTTGGGTGGACATCCGCTTCATCTCCCTCAAAGATCAAGGATGACTCATGATCTCTGAGGCCATAGTCACATGTGGTGGT

16►aProAsnLeuLeuGl yGl yProSer ValPhellePheProProLysI eLysAspValLeuMetI eSerLeuSerProIleValThrCysValVal

PvuII (835)

801 GATGTGAGCGAGGTGACCGATGTCAGCTCAGCTGGTTGTGAACACAGTGGAAAGTACACACAGCTCAGACACAAACCCATAGAGGGATTACAACA

50►AspValSerGl uAspAspProAspValGl nIeSerTrpPheValAsnAsnValGl uValHisThrAl aGl nThrGl nThrHi sArgGl uAspTyrAsnS

901 GTACTCTCGGGTGTCAAGTGCCTCCCTCCACGACCAAGGACTGGATGAGTGGCAAGGATTCAATGCAAGGTCACAACAAAGACCTCCAGGCC

83►eThrLeuArgValValSerAl aLeuProl IeGl nHi sGl nAspTrpMetSerGl yLysGl uPheLysCysLysValAsnAsnLysAspLeuP roAl aPr

1001 CATCGAGAGAACCATCTAAACCCAAAGGGTCAAGAGCTCCACAGGTATATGCTTGCCTCACCAGAAGAAGAGATGACTAAGAACACGGTACT

116►ol IeGl uArgThr I eSerLysProLysGl ySerValArgAl aProGl nValTyrValLeuProProGl uGl uGl uMetThrLysLysGl nValThr

1101 CTGACCTGCATGGTCACAGACTTCATGCCAGAACGACTTACGTTAGGGAGTGGACCAACAACGGAAAACAGAGCTAAACTACAAGAACACTGAACAGTCC

150►LeuThrCysMetValThrAspPheMetProGl uAspI IeTyrValGl uTrpThrAsnAsnGl yLysThrGl uLeuAsnTyrLysAsnThrGl uProValI

BsrGI (1225)

1201 TGGACTCTGATGGTCTTACTTCATGACAGACTGAGAGTGGAAAAGAAGAACCTGGTGGAAAGAAATAGCTACTCTGTTAGTGGTACAGGAGGG

183►euAspSerAspGl ySerTyrPheMetTyrSerLysLeuArgValGl uLysLysAsnTrpValGl uArgAsnSerTyrSerCysSerValValHi sGl uGl

MscI (1366)
NheI (1358)

1301 TCTGCACAATCACACACGACTAACAGCTTCTCCGGACTCCGGTAAATGAGCTAGCTCAGCTGGCAGACATGATAAGATACATTGATGAGTTGGAC

216►yLeuHi sAsnHi sHi sThrThrLysSerPheSerArgThrProGl yLys***

HpaI (1498)

1401 AAACACAACTAGAATGCAGTAAAAATGCTTTAGGTGAAATTGTGATGCTATTGTTATTGTAACCATTATAAGCTGCAATAAACAGTTAA

AsI (1595)
XmnI (1596)

1501 CAACAACAATTGATTCTATTTATGTTTCAAGTTCAAGGGGAGGTGGAGGTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGATTAAATT

1601 CTTAAATACAGCATAGCAAACCTTAACCTCAAATCAAGCTACTTGAATCTTCTGAGGGATGAATAAGGCATAGGCATCAGGGCTGTTGCC

1701 ATGTGCATTAGCTTTGAGCCTCACCTCTTCAAGGGTTAAGATATAGTGTATTGCCAGGTTGAACAGCTCTCATTTCTATGTTTA

SwaI (1852)

1801 AATGCACTGACCTCCACATCCCTTTAGTAAATATTCAAGAAATAATTAAACATCATTGCAATGAAATAATGTTTATTAGGCAGAACCT

1901 AGATGCTCAAGGCCCTCATATAATCCCCAGTTAGTAGTGGACTAGGGACAAAGAACCTTAATAGAAATTGGACAGCAAGAACGGCT

2001 AGCTTATCCTCAGTCTGCTCTGCCAACAAAGTGCAGCGACTGGCGGGGGCTCGCAGGGCGAACCTGGCTGCTGCCGATCTCG

125►•••AspGl nGl uGl uAl aVal PheHi sVal CysAsnGl yAl aProAspArgLeuAl aPheGl uArgGl yTrpProGl nGl uGl yI eGl uT

2101 GTCATGGCCGGCCGGAGGCCTCCGGAAAGTCTGAGCACGACCTCCGACACTCGCGTACAGCTCGTCCAGGGCGCACCCACCCAGGCCAGG

94►hr MetAl aProGl ySerAl aAspArgPheAsnThrSerValValGl uSerTrpGl uAl aTyrLeuGl uAspLeuGl yArgVal TrpValTrpAl aleuTh

XmaI (2291)
SmaI (2293)

2201 TGTGTCGGCACCACTGGCTTGGACCGCGCTGATGAACAGGGTCAGCTCGCCGGACACAGGGCGAACGTCTCCACGAAGTCCGGAGAA

61►rAsnAspProValValGl nAspGl nValAl aSerI IePheLeuThrValAspAspArgValValGl yAl aPheAspAspGl uValPheAspArgSerPhe

2301 CCCGAGCCGGTGGCTCCAGAACCTGACCGCTCCGGCGACGTCGCGCGGGTGGACACGGAACGGCACTGGCAACTTGGCCATGATGGCTCCTCtgc

28►Gl yLeuArgAspThrTrpPheGl uValAl aGl yAl aValAspArgAl aThrLeuValProValAl aSerThrLeuLysAl aMet

AsI (2479) PstI (2501)

2401 agggagggaaagagaagaaggtagtacaattgCTATAGTGGAGTTATTATCATGTCAGATATACTATGCAATGATTAATTGCAAATAGGGCTG

HindIII (2605)

2501 AggttcatagtccactttctgcactgccccatctcctgcccacccttccaggcatagacagtcaacttgactacCAAAC

ACTCACAGGAGGGAGAAG

2601 GCAGAACGCTTGAGACAGACCCGGGGACCGCCGAUTGCGAGGGACGGCTAGGGCTTCTTATGGTGCCTGGGAGGGCCTGGAGGGCTC

←

StuI (2707)
Eco147I (2707)

2701 GGGGAGGCCAGGCCAATCTCGGTGGCAGGAGGCAGGGCCGAAGGCCGTGCCGACCAATCCGGAGCACATAGGAGTCTCAGCCCCCGCCCCAAAG

BspEI (2763)

2801 CAAGGGGAAGTCACGCCCTGTAGGCCAGCGTTGAAATGGGGCTGGGGGGTGGGCCCTGACTAGTC**C**AAAACAAACTCCCATTGACGTCAA

SpeI (2870)

2901 TGGGGTGGAGACTTGGAAATCCCGTGAGTC**A**AAACCGCTATCCACGCCATTGATGTACTGCCAAACCGCATCATGTAATAGCGATGACTAAC

SnaBI (3000)

3001 GTAGATGACTGCCAAGTAGGAAAGTCCCATAAGGTATGACTGGCATAATGCCAGGCCATTACCGTATTGACGTCAATAGGGCGTACTT

NdeI (3104)

3101 GGCATATGATACACTTGATGACTGCCAAGTGGCAGTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCATTGGCGTTACTATGGAA

PacI (3293)

3201 CATACGTATTATTGACGTCAATGGCGGGGTCGTTGGCGGTAGCCAGGCAGGCCATTACCGTAAGTTATGTAACGCCCTGCAGGTTAATTAGAAC

PstI (3286)
SdaI (3286)
BspLU1II (3299)

3301 ATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAGGCCGTTGCTGCCGTTTCCATAGGCTCGCCCCCTGACGAGCATCACAAATC

3401 GACGCTCAAGTCAGAGGTGGCAAACCCGACAGGACTATAAGATAACCAGCGTTCCCTGGAAAGCTCCCTGCGCTCTCTGTTCCGACCCGCC

3501 GCTTACCGGATACTTGTCCGCCCTTCTCCCTGGAAAGCGTGGCGTTCTCATAGCTCACGCTGTAGGTATCTAGTTGGTAGGTCTGTTCCGCTCC

3601 AAGCTGGCTGTGTCACGAACCCCCGTTAGCCGCTGCCCTATCCGTAACTATCGTCTGAGTCCAACCCGTAAGACACGACTATCGC

3701 CACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGAGCGGTGCTACAGAGTTCTGAAGTGGTGGCTAACTACGGTACACTAGAAG

3801 AACAGTATTGGTATCTGCGCTGCTGAAGCCAGTTACCTCGAAAAAGAGTTGGTAGCTTGATCCGCAACAAACACCACGCTGGTAGCGGTGGT

3901 TTTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAAGGATCTCAAGAAGATCCTTGATCTTCTACGGGCTGACGCTCAGTGGAACGAAACT

PacI (4033) **Swal (4041)** **NotI (4049)**

4001 CACGTTAAGGGATTGGTCATGGCTAGTTAATTAAACATTAAATCAGCGGCCAATAAAATATCTTATTTATTACATCTGTTGGTTTTTG

4101 TGTGAATCGTAACAAACATACGCTCTCCATCAAACAAACAAACGAAACAAACAAACTAGCAAATAGGCTGCCCCAGTGCAGGTGCCAGAACAT

4201 TTCTCTATCGAA