

pFUSE-mIgG2A-Fc1

Plasmid containing a mouse IgG2A Fc region

Catalog # pfuse-mg2afc1

For research use only

Version # 08F06-SV

PRODUCT INFORMATION

Content:

- 20 µg of pFUSE-mIgG2A-Fc1 plasmid provided as lyophilized DNA
- 4 pouches of *E. coli* Fast-Media® Zeo (2 TB and 2 Agar)

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 *E. coli* Fast-Media® Zeo at room temperature. Fast-Media® pouches are stable 18 months when stored properly.

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-Fc plasmids allow the secretion of Fc-Fusion proteins (when Fc portion is fused to a naturally secreted protein). 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.

In vivoGen 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.
- **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 *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.

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

Toll free (US): 888-457-5873
Outside US: (+1) 858-457-5873
E-mail: info@invivogen.com
Website: www.invivogen.com



3950 Sorrento Valley Blvd. Suite 100
San Diego, CA 92121 - USA

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.

Selection of bacteria with *E. coli* Fast-Media®

Fast-Media® is a **fast and convenient** way to prepare liquid and solid media for bacterial culture by using only a microwave. Fast-Media® is a TB (liquid) or LB (solid) based medium that already contains the antibiotic. Fast-Media® Zeo is available separately: #fas-zn-l (liquid), #fas-zn-s (agar).

- 1- Pour the contents of a Fast-Media® pouch into a clean borosilicate glass bottle or flask.
- 2- Add 200 ml of distilled water to the flask
- 3- Heat in a microwave on MEDIUM power setting (about 400Watts), until bubbles start appearing (approximately 3 minutes). **Do not heat a closed container. Do not autoclave Fast-Media®.**
- 4- Swirl gently to mix the preparation. **Be careful, the bottle and media are hot, use heatproof pads or gloves and care when handling.**
- 5- Reheat the media for 30 seconds and gently swirl again. Repeat as necessary to completely dissolve the powder into solution. But be careful to avoid overboiling and volume loss.
- 6- Let agar medium cool to 45°C before pouring plates. Let liquid media cool to 37°C before seeding bacteria.

Note: Do not reheat solidified Fast-Media® as the antibiotic will be permanently destroyed by the procedure.

RELATED PRODUCTS

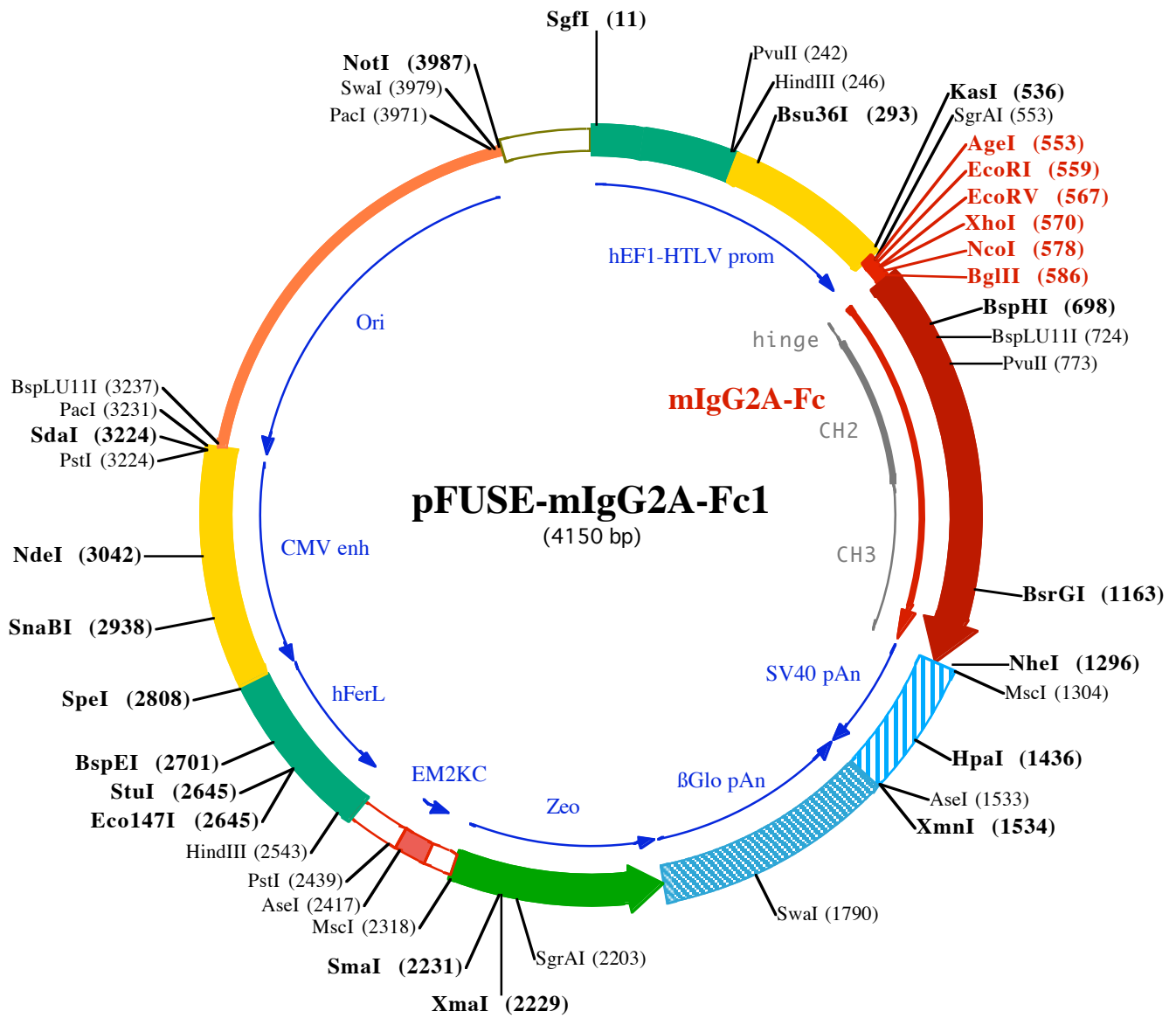
Product	Catalog Code
Zeocin™	ant-zn-1
Fast-Media® Zeo TB	fas-zn-l
Fast-Media® Zeo Agar	fas-zn-s

TECHNICAL SUPPORT

Toll free (US): 888-457-5873
Outside US: (+1) 858-457-5873
E-mail: info@invivogen.com
Website: www.invivogen.com



3950 Sorrento Valley Blvd. Suite 100
San Diego, CA 92121 - USA



SgfI (11)
1 GGATCTGCGATCGTCCCGTCCCGTCCAGTGGGCGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGTGGCAATTGAACGGTGCCTA
101 GAGAAGGTGGCGGGGTAACCTGGGAAAGTGATGTCTGTACTGGCTCCGCTTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGGC

HindIII (246)
PvuII (242)
Bsu36I (293)
201 GTGAACGTTCTTTTTCGCAACGGGTTTGGCCGAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACGCGCCCGCCCTACCTGAGGCC
301 GCCATCCACGCGGGTTGAGTGCCTCTGCCGCTCCCGCTGTGGTGCCTCTGAACCTGCTCCGCGTCTAGGTAAGTTAAAGCTCAGGTCGAGACC
401 GGGCTTTGTCCGGCGCTCCCTGGAGCTACCTAGACTCAGCGGCTCTCCACGCTTGGCTGACCTGCTTGTCTCAACTCTACGCTTTGTTCGTTT

KasI (536)
AgeI (553)
EcoRI (559)
XhoI (570)
BglII (586)
SgrAI (553)
EcoRV (567)
NcoI (578)
501 TCTGTTCTGGCGGTTACAGATCCAAGCTGTGACCGGCGCTACCTGAGATCACCGGTGAATTCGATATCTCGAGCACCATGGTTAGACTCTCCAGAGGG
1►ProArgGIy

BspHI (698)
601 CCCACAATCAAGCCTGTCTCCATGCAATGCCAGCACCTAACCTTTGGGTGGACCTCCGCTTTCATCTTCCCTCAAAGATCAAGGATGTACTCA
4►ProThrIleLysProCysProP roCysLysCysProAl aProAsnLeuLeuGI yGI yProSerValPheIlePheProProLysIleLysAspValLeuM

BspLU111 (724)
PvuII (773)
701 TGATCTCCCTGAGCCCATAGTCACATGTGTGGTGGTGGATGTGAGCGAGGATGACCCAGATGTCCAGATCAGCTGGTTTGTGAACAACGTGGAAGTACA
37►eIleSerLeuSerProIleValThrCysValValValAspValSerGIuAspAspProAspValGlnIleSerTrpPheValAsnAsnValGluValHi
801 CACAGCTCAGACAAAACCCATAGAGAGGATTACAACAGTACTCTCCGGTGGTGCAGTGCCTCCCATCCAGCACCAGGACTGGATGAGTGGCAAGGAG
70►sThrAlaGlnThrGlnThrHisArgGIuAspTyrAsnSerThrLeuArgValValSerAlaLeuProIleGlnHisGlnAspTrpMetSerGIyLysGIu
901 TTCAAATGCAAGGTCAACAACAAGACCTCCAGCGCCATCGAGAGAACCATCTCAAAACCAAGGGTGCAGTAAGAGCTCCACAGGTATATGTCTTGC
104►PheLysCysLysValAsnAsnLysAspLeuProAlaProIleGIuArgThrIleSerLysProLysGIySerValArgAlaProGlnValTyrValLeuP
1001 CTCCACAGAAGAAGAGATGACTAAGAAACAGTCACTGACCTGCATGGTCACAGACTTCATGCCTGAAGACATTTACGTGGAGTGGACCAACAACGG
137►roProProGIuGIuGIuMeThrLysLysGIuValThrLeuThrCysMetValThrAspPheMetProGIuAspIleTyrValGIuTrpThrAsnAsnGI

BsrGI (1163)
1101 GAAAACAGAGCTAACTACAAGAACTGAACAGTCTGGACTCTGATGGTTCTTACTTCATGTACAGCAAGCTGAGAGTGGAAAAGAAGAACTGGGTG
170►yLysThrGIuLeuAsnTyrLysAsnThrGIuProValIleuAspSerAspGIySerTyrPheMetTyrSerLysLeuArgValGIuLysLysAsnTrpVal

NheI (1296)
1201 GAAAGAAATAGCTACTCTGTTCAGTGGTCCAGAGGCTGCACAATCACACAGACTAAGAGCTTCTCCCGACTCCGGTAATGAGCTCAGCTAG
204►GIuArgAsnSerTyrSerCysSerValValHisGIuGIyLeuHisAsnHisHisThrThrLysSerPheSerArgThrProGIyLys●●●

MscI (1304)
1301 CTGGCAGACATGATAAGATACATTGATGAGTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTGTGAAATTTGTGATGCTATTGCT

HpaI (1436)
1401 TTATTTGAACATTATAAGCTGCAATAAACAAGTTAAACAACAACATTCATTCATTTTATGTTTCAGGTTCCAGGGGAGGTGGGAGGTTTTTAA

AseI (1533)
XmnI (1534)
1501 GCAAGTAAACCTCTACAAATGTGGTATGGAATTAATCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTG
1601 AGGGATGAATAAGGCATAGGCATCAGGGCTGTGCCAATGTGCATTAGCTGTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTTCT

Swal (1790)
1701 CCAAGTTTGAAGTACTCTTCAATTTCTTATGTTTTAAATGACTGACCTCCACATTCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCA
1801 TTGCAATGAAAATAAATGTTTTTATTAGCGAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAGGAA
1901 CCTTTAATAGAAATGGACAGCAAGAAGCGAGCTTCTAGCTTATCTCAGTCTGCTCTGCCACAAAGTGCACGAGTTGCCGCGCGGGTCCGCGCA
125►●●AspGlnGIuGIuAlaValPheHisValCysAsnGIyAlaProAspArgLe
2001 GGGCAACTCCGCCCCACGGCTGCTCGCCGATCTCGGTATGGCGGCCGGAGGCGTCCCGAAAGTTCGTGGACACGACTCCGACACTCGGCGTA
107►AlaPheGIuArgGIyTrpProGlnGIuGIyIleGIuThrMetAlaProGIySerAlaAspArgPheAsnThrSerValValGIuSerTrpGIuAlaTyr
2101 CAGCTCGTCCAGGCGCGCACCCACACCCAGGCCAGGCTGTGTCCGGCACCTGGTCTGGACCGCGCTGATGAACAGGGTCACTGCTCCCGACC
74►LeuGIuAspLeuGIyArgValTrpValTrpAlaLeuThrAsnAspProValValGlnAspGlnValAlaSerIlePheLeuThrValAspAspArgValV

XmaI (2229)
SmaI (2231)
2201 ACACCGCGAAGTCTCTCCACGAAGTCCCGGAGAACCCGAGCGGTGCGTCCAGAATCGACCGCTCCGCGCAGCTCGCGCGCGGTGAGCACCGGAA
40►AlaGIyAlaPheAspAspGIuValPheAspArgSerPheGIyLeuArgAspThrTrpPheGIuValAlaGIyAlaValAspArgAlaThrLeuValProVa
2301 CGGCACTGGTCAACTGGCCATGATGGCTCTCctgtcaggagaggaagagaagaaggttagtacaattgCTATAGTGAGTTGTATTACTATGCGAGA
7►AlaIaSerThrLeuLysAlaMe t
2401 TATACTATGCCAATGATTAATTGTCAAACCTAGGGCTGCAgggttcatagtgcacttttctgcactgccccatctcctgccccctttccaggcata

HindIII (2543)
2501 gacagtcagtgacttacCAAACCTACAGGAGGGAGAAGCGAAGCTTGAGACAGACCCCGGGACCGCCAAGTGCAGGGGACGTGGCTAGGGCGGT

StuI (2645)
Eco147I (2645)
2601 TCTTTTATGTTGCGCCGCCCTCGAGGCGAGGCGCTCGGGGAGGCTAGCGGCCAATCTGCGGTGGCAGGAGCGGGCCGAAGGCGGTGCTGACCAA

2701 **BspEI (2701)**
TCCGGAGCACATAGGAGTCTCAGCCCCCGCCCCAAAGCAAGGGGAAGTCACGCGCCTGTAGGCCAGCGTGTGTGAAATGGGGCTTGGGGGGTTGG

2801 **SpeI (2808)**
GGCCCTGACTAGTCAAACAACAACTCCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTA CTGC

2901 **SnaBI (2938)**
CAAAACCCGATCATCATGGTAATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCATAAGGTCATGTA CTACTGGGCATAATGCCAGGCGG

3001 **NdeI (3042)**
GCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGCAGTTTACCGTAAATACTCCACCCATTG

3101 ACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGCGGTCAGCCAGGCGGGCCATTT

3201 **SdaI (3224)** Pacl (3231) PstI (3224) BspLU11I (3237)
ACCGTAAGTTATGTAACGCCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTT

3301 CCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCCT

3401 GGAAGCTCCCTCGTGGCTCTCCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCAC

3501 GCTGTAGGTATCTCAGTTCGGGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCCAGCCGACCGCTGCCTTATCCGGTAAC TA

3601 TCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAG

3701 TTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCT

3801 CTTGATCCGGCAAACAACCACCGCTGGTAGCGGTGGTTTTTTTGTGCAAGCAGCAGATTACGGCAGAAAAAAGGATCTCAAGAAGATCCTTTGAT

3901 CTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAA

4001 TATCTTTATTTTCATTACATCTGTGTGGTTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACAACTAGCAAA

4101 ATAGGCTGTCCCCAGTGAAGTGCAGGTGCCAGAACATTTCTCTATCGAA