

# pFUSE-hIgG4-Fc2

Plasmid designed for the construction of Fc-Fusion proteins

Catalog # pfuse-hg40fc2

For research use only

Version 20K05-MM

## PRODUCT INFORMATION

### Content:

- 20 µg of pFUSE-hIgG4-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. In humans, there are four isotypes: IgG1, IgG2, IgG3 and IgG4. The Fc region mediates effector functions, such as antibody-dependent cellular cytotoxicity (ADCC) and complement-dependent cytotoxicity (CDC). IgG isoforms exert different levels of effector functions increasing in the order of IgG4<IgG2<IgG1<IgG3.

## PLASMID FEATURES

- **hIgG4-Fc (human):** 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. Human IgG4 displays low ADCC and no CDC, and therefore is the most suitable for diagnostic imaging or blocking molecular interactions<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*.
- **BGlo pAn:** The human beta-globin 3'UTR and polyadenylation sequence allows efficient arrest of the transgene transcription<sup>5</sup>.

1. Kim SJ. et al., 2005. Antibody engineering for the development of therapeutic antibodies. *Mol. Cells*. 20(1):17-29.
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

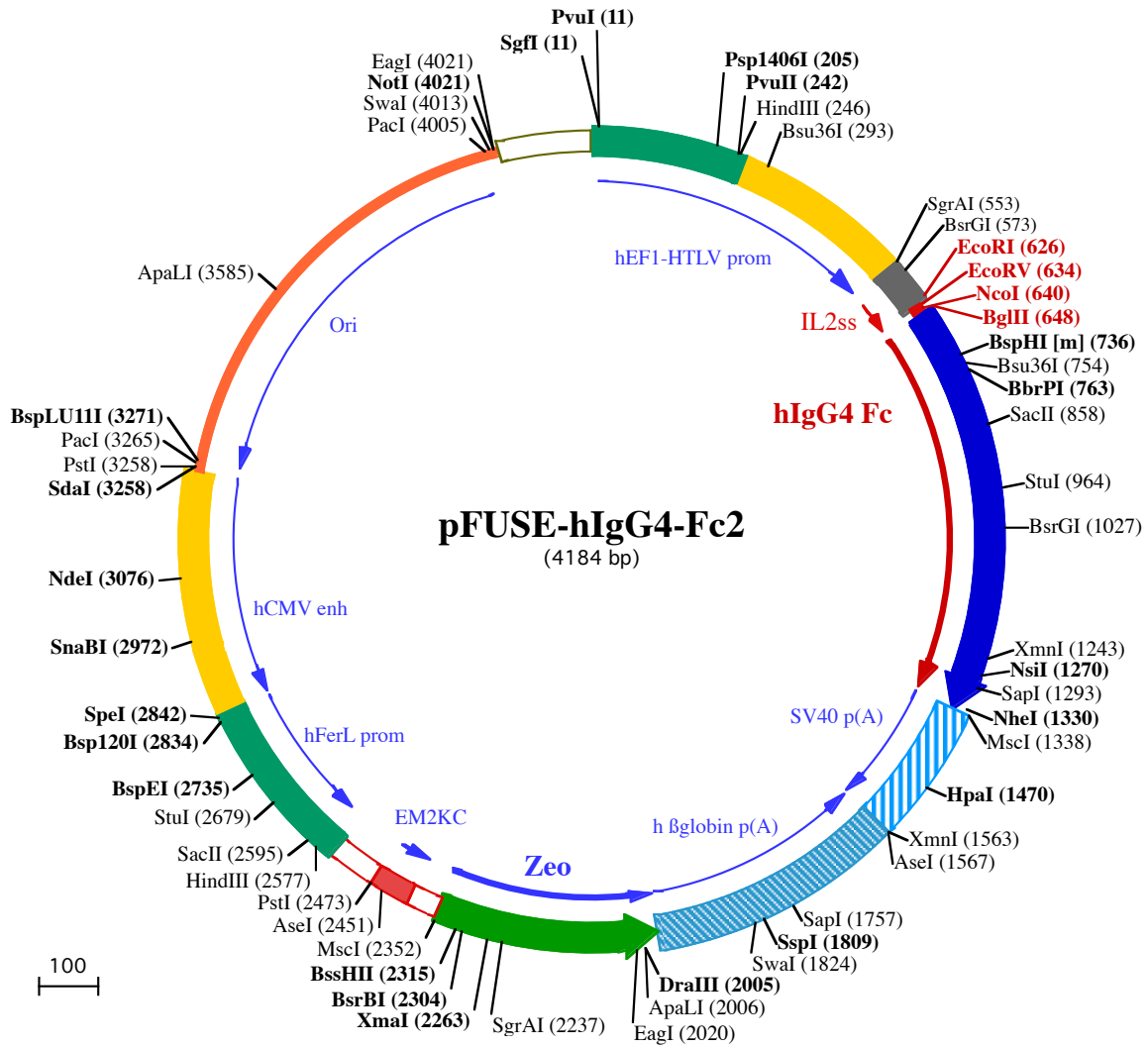
InvivoGen USA (Toll-Free): 888-457-5873

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)



**PvuI (11)**  
**SgfI (11)**  
 1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA  
 101 GAGAAAGTGGCGCGGGTAAACTGGAAAAGTATGTCGTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

---

**HindIII (246)**  
**Psp1406I (205)** **PvuII (242)** **Bsu36I (293)**  
 201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACCGCGCCGCCCTACCTGAGGCC  
 301 GCCATCCACGCCGGTTGAGTCGCGTTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAATGCGTCCCGCTTAGGTAAGTTTAAAGCTCAGGTCGAGACC  
 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTCCTGACCTGCTTGTCTAACTCTACGCTTTTGTTCGTTT

---

**SgrAI (553)** **BsrGI (573)**  
 501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGCGAAGGAGGGCCACCATGTACAGGATGCAACTCCTGTCTTGCA  
 1▶ M Y R M Q L L S C

---

**EcoRV (634)**  
**EcoRI (626)** **NcoI (640)** **BglIII (648)**  
 601 TTGCACTAAGTCTTGCACTTGTACGAATTCGATATCGGCCATGGTTAGATCTCCCCATGCCATCATGCCAGCACCTGAGTTCCTGGGGGACCATC  
 10▶ I A L S L A L V T N S 1▶ P P C P S C P A P E F L G G P S

---

**BbrPI (763)**  
**BspHI [m] (736)** **Bsu36I (754)**  
 701 AGTCTTCTGTTCCTCCCAAAACCAAGGACTCTCATGATCTCCGGACCCCTGAGGTACGTCAGTGGTGGTGGACGTGAGCCAGGAAGACCCCGAG  
 16▶ V F L F P P K P K D T L M I S R T P E V T C V V V D V S Q E D P E

---

**SacII (858)**  
 801 GTCCAGTTCAACTGGTACGTGGATGGCGTGGAGGTGCATAATGCCAAGACAAAGCCGCGGGAGGAGCAGTTCAACAGCACGTACCGTGTGGTACGCTCC  
 50▶ V Q F N W Y V D G V E V H N A K T K P R E E Q F N S T Y R V V S V

---

**StuI (964)**  
 901 TCACCGTCTGCACCAGGACTGGCTGAACGGCAAGGAGTACAAGTGAAGGTCTCAACAAAGGCCTCCCGTCTCCATCGAGAAAACCATCTCCAAAGC  
 83▶ L T V L H Q D W L N G K E Y K C K V S N K G L P S S I E K T I S K A

---

**BsrGI (1027)**  
 1001 CAAAGGGCAGCCCGAGAGCCACAGGTGTACACCTGCCCCATCCCAGGAGGAGATGACCAAGAACCAGGTGAGCCTGACCTGCCTGGTCAAAGGCTTC  
 116▶ K G Q P R E P Q V Y T L P P S Q E E M T K N Q V S L T C L V K G F  
 1101 TACCCAGCGACATCGCCGTGGAGTGGGAGAGCAATGGGAGCCGGAGACAACACTACAAGACCACGCCTCCCGTGTGGACTCCGACGGCTCCTTCTTCC  
 150▶ Y P S D I A V E W E S N G Q P E N N Y K T T P P V L D S D G S F F

---

**XmnI (1243)** **NsiI (1270)** **SapI (1293)**  
 1201 TCTACAGCAGGCTAACCTGGACAAGAGCAGGTGGCAGGAGGGGAATGCTTCTCATGCTCCGTGATGCATGAGGCTCTGCACAACCACTACACACAGAA  
 183▶ L Y S R L T V D K S R W Q E G N V F S C S V M H E A L H N H Y T Q K

---

**MscI (1338)**  
**NheI (1330)**  
 1301 GAGCCTCTCCCTGTCTCIGGGTAAATAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAA  
 216▶ S L S L S L G K •

---

**HpaI (1470)**  
 1401 ATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGTGAATAAACAAGTTAACAACAACAAATTGCATTATTTTATGTTT

---

**AseI (1567)** **XmnI (1563)**  
 1501 CAGGTTACAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATTAATTCTAAAATACAGCATAGCAAACTTTAAC  
 1601 CTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCAC

---

**SapI (1757)**  
 1701 TTCTTTCATGGAGTTTAAGATATAGTGTATTTTCCCAAGGTTTGAAGTACTGCTCTTCTTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTT

---

**SspI (1809)** **SwaI (1824)**  
 1801 TAGTAAAAATATTCAGAAATAAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCC

---

**DraIII (2005)**  
 1901 CCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTAAATAGAAATTTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTCTGCC  
 125▶ • D Q E E A

---

**ApaLI (2006)** **EagI (2020)**  
 2001 ACAAAGTGCAGCAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTCATGGCCGGCCCGAGGCGTCCCGGA  
 118▶ V F H V C N G A P D R L A F E R G W P Q E G I E T M A P G S A D R F  
 2101 AGTTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCGCGCACCCACACCCAGGCCAGGGTGTGTCGGCACCACCTGGTCTGGAC  
 85▶ N T S V V E S W E A Y L E D L G R V W V W A L T N D P V V Q D Q V

---

**SgrAI (2237)** **XmaI (2263)** **BsrBI (2304)**  
 2201 CGCGCTGATGAACAGGGTCACGTCGTCGGGACACACCGCGGAAGTCGTCTCCACGAAGTCCCGGAGAACCCGAGCCGGTCCGTCAGAACTCGACC  
 52▶ A S I F L T V D D R V V G A F D D E V F D R S F G L R D T W F E V

2301 **BssHII (2315)** **MscI (2352)**  
 18 GCTCCGGCGACGTCGCGCGCGGTGAGCACCGGAACGGCACTGGTCAACTTGGCCATGATGGCTCCTCctgtcaggagaggaagagaagaaggtagtac  
 18 A G A V D R A T L V P V A S T L K A M

2401 **AseI (2451)** **PstI (2473)**  
 aattgCTATAGTGAGTTGTATTATACTATGCAGATATACTATGCCAATGATTAATTGTCAAAC TAGGGCTGC Agggttcatagtgccacttttctgcac

2501 **HindIII (2577)** **SacII (2595)**  
 tgcctcatctcctgccaccctttccaggcatagacagtcagtgacttacCAAAC TACAGGAGGAGAAGGCAGAAAGCTTGAGACAGACCCGCGGGAC

2601 **StuI (2679)**  
 CGCCGAAC TCGAGGGGACGTGGCTAGGGCGGCTTCTTTTATGGTGC CGCCCTCGGAGGCAGGGCGCTCGGGGAGGCCTAGCGGCCAATCTGCGGTG

2701 **BspEI (2735)**  
 GCAGGAGCGGGGCCGAAGGCCGTGCCTGACCAATCCGGAGCACATAGGAGTCTCAGCCCCCGCCCAAAGCAAGGGGAAGTCACGCGCCTGTAGCGCC

2801 **SpeI (2842)** **Bsp120I (2834)**  
 AGCGTGTGTGAAATGGGGGCTTGGGGGGTTGGGGCCCTGACTAGTCAAACCAAAC TCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGA

2901 **SnaBI (2972)**  
 GTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCC

3001 **NdeI (3076)**  
 CATAAGGTCATGTACTGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATAATGATACACTTGATGTACTGCCA

3101 **NdeI (3076)**  
 AGTGGGCAGTTTACCCTAAATACTCCACCCATTGACGTCAATGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCAATATTGACGTCAATGGGGC

3201 **PacI (3265)** **PstI (3258)** **SdaI (3258)** **BspLU11I (3271)**  
 GGGTTCGTTGGCGGTACAGCCAGGCGGGCCATTACCCTAAGTTATGTAACGCCTGCAGGTTAATTAAGAACATGTGAGCAAAAAGCCAGCAAAAAGGCCA

3301 **BspLU11I (3271)**  
 GGAACCGTAAAAAGCCGCTTGTGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCC

3401 **BspLU11I (3271)**  
 GACAGGACTATAAAGATACCAGGCGTTCCCCCTGGAAGCTCCCTCGTGCGCTCTCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTC

3501 **ApaLI (3585)**  
 CCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGATAGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGC

3601 **ApaLI (3585)**  
 TTCAGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGAT

3701 **ApaLI (3585)**  
 TAGCAGAGCGAGGTATGTAGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTG

3801 **ApaLI (3585)**  
 AAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGCAAGCAGCAGATTACGC

3901 **ApaLI (3585)**  
 GCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAAACGAAAAC TACGTTAAGGGATTTTGGTCATGGCTAG

4001 **EagI (4021)** **PacI (4005)** **SwaI (4013)** **NotI (4021)**  
 TTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTTCATTACATCTGTGTGTTGGTTTTTTTGTGTAATCGTAACTAACATACGCTCTCC

4101 **NotI (4021)**  
 ATCAAAACAAAACGAAACAAAACAAACTAGCAAATAGGCTGTCCCCAGTGAAGTGCAGGTGCCAGAACATTTCTCTATCGAA