

# pFUSE-hIgG1e11-Fc1

Plasmid containing a human engineered IgG1 Fc region

Catalog # pfc1-hg1e11

For research use only

Version 20K05-MM

## PRODUCT INFORMATION

### Content:

- 20 µg of pFUSE-hIgG1e11-Fc1 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-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 protein A 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. Human IgG1 displays high ADCC and CDC, and is the most suitable for therapeutic use against pathogens and cancer cells.

Under certain circumstances, for example when depletion of the target cell is undesirable, abrogating effector functions is required. On the contrary, in the case of antibodies intended for oncology use, increasing effector functions may improve their therapeutic activity<sup>1</sup>. Modifying effector functions can be achieved by engineering the Fc regions to either improve or reduce their binding to FcγRs or the complement factors. Amino acid substitutions have been made in the human IgG1 Fc region in order to increase or reduce its ADCC and CDC.

## PLASMID FEATURES

- **hIgG1e11-Fc (human IgG1 engineered Fc):** 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 binds to neonatal FcR (FcRn), a receptor expressed on the surface of endothelial cells. This interaction, which is pH-dependent, protects the IgG from lysosomal degradation thus mediating the serum persistence of IgG antibodies. The human IgG1 Fc domain was engineered by introducing mutations in the FcRn binding sites leading to higher FcRn binding affinity at pH 6.0<sup>2</sup>. The engineered pFUSE-hIgG1e11-Fc1 contains three amino acid substitutions: M252Y, S254T and T256E.
- **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1α (EF-1α) core promoter<sup>3</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>4</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.
- **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>5</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>6</sup>.

1. Carter PJ., 2006. Potent antibody therapeutics by design. Nature Reviews Immunology. Nat Rev Immunol. 6, 343-357.
2. Oganessian V. *et al.*, 2009. Structural characterization of a human Fc fragment engineered for extended serum half-life. Molec Immunol 46: 1750-1755.
3. Kim DW *et al.* 1990. Use of the human elongation factor 1 alpha promoter as a versatile and efficient expression system. Gene 91(2):217-23.
4. 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.
5. 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.
6. 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](#)

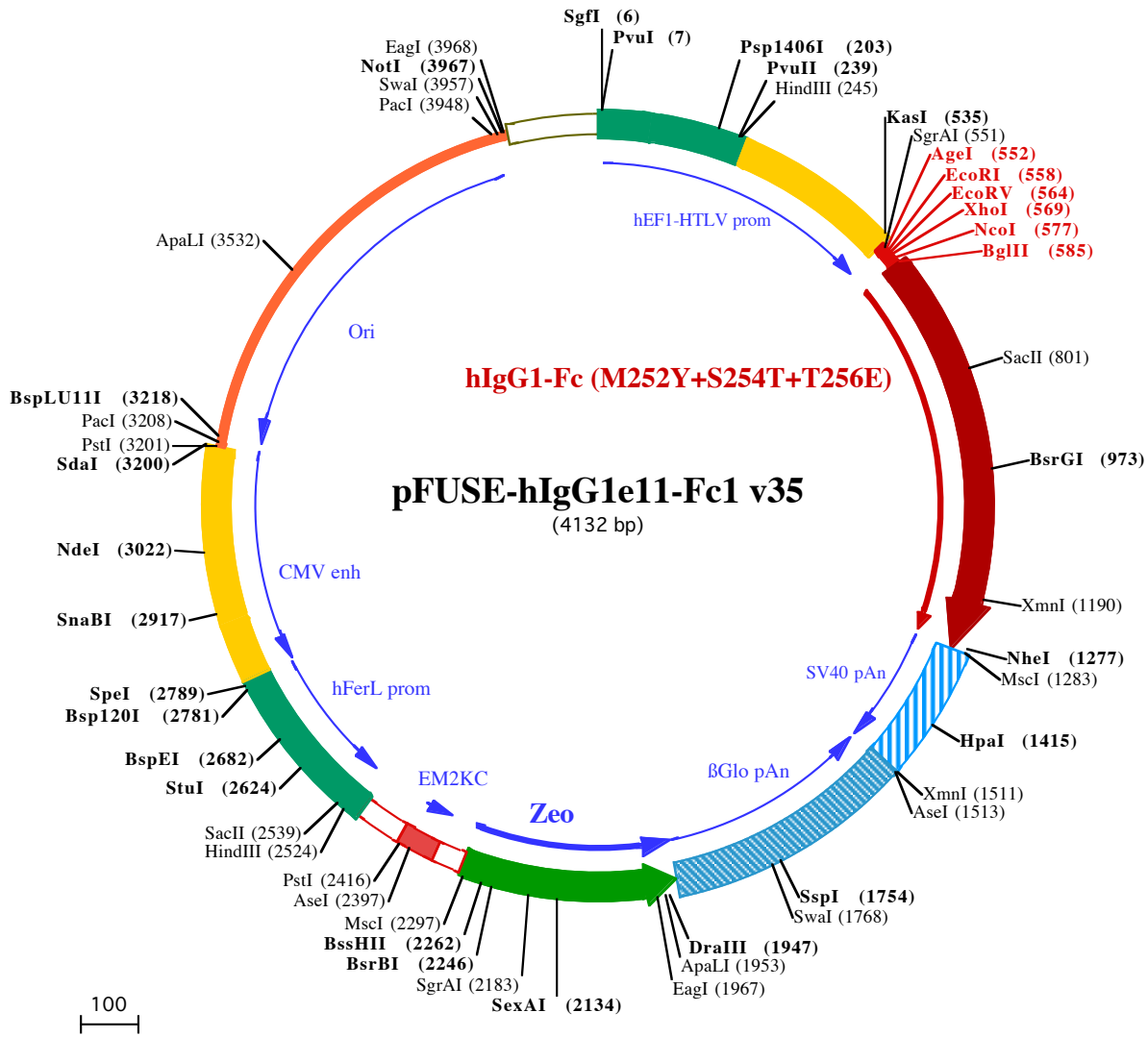
InvivoGen USA (Toll-Free): 888-457-5873

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 (7)  
SgfI (6)  
1 GGATCTGCGATCGCTCCGGTCCCGTCCAGTGGGACAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGTGGCAATTGAACGGTGCCTA  
101 GAGAAAGTGGCGGGGTAAGTGGGAAAGTGATGCTGTACTGGCTCCGCCCTTTTCCGAGGGTGGGGGAGAACGTATATAAGTGCAGTAGTCGCC

HindIII (245)  
Psp1406I (203) PvuII (239)  
201 GTGAACGTTCTTTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTCGAGGGGTCGCATCTCTCTTCACGCGCCCGCCCTACCTGAGGCC  
301 GCCATCCACGCCGGTTGAGTCGCTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAATGCGTCCGCGCTAGGTAAGTTTAAAGCTCAGTGCAGAGCC  
401 GGGCCTTTGTCCGGCGCTCCCTTGAGGCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

EcoRI (558) AgeI (552) XhoI (569) BglII (585)  
KasI (535) SgrAI (551) EcoRV (564) NcoI (577)  
501 TCTGTTCTGCCCGTTACAGATCCAAGCTGTGACCGGCGCCTACCTGAGATCACCGGTAATTCGATATCTCGAGCACCATGGTTAGATCTGACAAAAC  
601 CACACATGCCACCGTGGCCAGCCTGAACCTCTGGGGGACCGTCACTCTTCTCTTCCCCAAAACCAAGGACACCTCTACATCAACCCGGGAA  
700 CCTGAGGTACATGCGTGGTGGTGGAGCTGAGCCAGAACCCCTGAGTCAAGTTCAACTGGTACGTGGACGGCGTGGAGGTGCATAATGCCAAGACAA  
37▶ P E V T C V V V D V S H E D P E V K F N W Y V D G V E V H N A K T  
SacII (801)  
800 AGCCGCGGGAGGAGCAGTACAACAGCAGCTACCGTGTGGTCAAGCTCCTCACCGTCTGCACCAGGACTGGCTGAATGGCAAGGAGTACAAGTCAAGGT  
70▶ K P R E E Q Y N S T Y R V V S V L T V L H Q D W L N G K E Y K C K V  
BsrGI (973)  
900 CTCCAACAAAGCCCTCCAGCCCCATCGAAGAACCTCTCAAAGCCAAAGGGCAGCCCGAGAACCACAGGTGTACACCTGCCCCATCCCGGGAG  
103▶ S N K A L P A P I E K T I S K A K G Q P R E P Q V Y T L P P S R E  
1000 GAGATGACCAAGAACCGGTGAGCTGACCTGGTCAAAGCTTATCCAGCGACATCGCCGTGGAGTGGAGAGCAATGGCGAGCGGAGAACCA  
137▶ E M T K N Q V S L T C L V K G F Y P S D I A V E W E S N G Q P E N  
XmnI (1190)  
1100 ACTACAAGACCAGCCTCCGTGTGGACTCCGACGGCTCTTCTTCTCTACAGCAAGCTCACCGTGGACAAGAGCAGGTGGCAGCAGGGGAACGTCTT  
170▶ N Y K T T P P V L D S D G S F F L Y S K L T V D K S R W Q Q G N V F  
MscI (1283)  
NheI (1277)  
1200 CTCATGCTCCGTGATGCACGAGGCTCTGCACAACACTACAGCAGAAGAGCCTCTCCCTGTCTCCGGGTAATGAGTGTAGCTGGCCAGACATGATAA  
203▶ S C S V M H E A L H N H Y T Q K S L S L S P G K •  
1300 GATACATTGATGAGTTTGGACAAACCACTAGAATGCAGTGAAGAAATGCTTTATTTGTGAATTTGTGATGCTATTGCTTTATTTGTAACATTAT

HpaI (1415)  
1400 AAGCTGCAATAAACAAGTTAAACAACAATTGCATTATTTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAACCTCTAC

AseI (1513) XmnI (1511)  
1500 AAATGTGGTATGGAATTAATTCTAAAAACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCAT  
1600 AGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTCTTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGGTTTGAAGTAC

SspI (1754) SwaI (1768)  
1700 TCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTATGATAAATATTAGAAATATTAGAAATAATTTAAATACATCATTGCAATGAAAAATAAT  
1800 GTTTTTTATTAGGCAGAAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAATTGG

ApaLI (1953) DraIII (1947) EagI (1967)  
1900 ACAGCAAGAAAGCGAGCTTCTAGCTTATCTCAGTCTGCTCCTCTGCCACAAAGTGCACGAGTTGCCGGCCGGGTGCGCGAGGGCGAACTCCCGCCCC  
2000 CACGGCTGCTCGCGATCTCGTTCATGGCCGGCCGGAGCGTCCCGAAGTTCGTGGACACGACCTCCGACCACTCGCGTACAGCTCGTCCAGCCGCG  
101▶ W P Q E G I E T M A P G S A D R F N T S V V E S W E A Y L E D L G R  
SexAI (2134) SgrAI (2183)  
2100 GCACCCACACCCAGGCCAGGTGTTGTCCGGACCACTGGTCTGGACCGCGTGATGAACAGGGTACGTCGTCGCCGACACACCGCGAAGTCTGTC  
68▶ V W V W A L T N D P V V Q D Q V A S I F L T V D D R V V G A F D D  
BsrBI (2246) BssHII (2262) MscI (2297)  
2200 CTCCACGAAGTCCCGGAGAACCCGAGCCGGTCCGAGAACCTGACCGCTCCGGCGACGTCGCGCGGTTGAGCACCAGGACCGCACTGGTCAACTTG  
35▶ E V F D R S F G L R D T W F E V A G A V D R A T L V P V A S T L K  
AseI (2397)  
2300 GCCATGATGGCTCCTcctgtcaggagaggaagagaagaagggttagtacaattgCTATAGTGTATTATACTATGCAGATATACTATGCCAATGAT  
1▶ A M

PstI (2416)  
2400 TAATTGTCAAAGTGGCTGCAggggttatagtgccacttttctgctactgccccatctcctgccccaccctttccaggcatagacagtcagtgacttac

HindIII (2524) SacII (2539)  
2500 CAAACTCACAGGAGGAGAAGGCAGAAGCTTGAGACAGACCCGCGGACCCGCAACTGCGAGGGGACGTGGCTAGGCGGCTCTTTTATGGTGCGCCG

StuI (2624) BspEI (2682)  
2600 GCCCTCGGAGGACGGGCTCGGGGAGGCTAGCGGCAATCTGCGGTGGCAGGAGCGGGGCCAAGGCCGTGCTGACCAATCCGGACACATAGGAG

SpeI (2789) Bsp120I (2781)  
2700 TCTCAGCCCCCGCCAAAGCAAGGGAAGTACCGCCCTGTAGCGCCAGCTGTTGTGAAATGGGGGTTGGGGGTTGGGCCCTGACTAGTCAAA  
2800 ACAAACTCCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCAT

SnaBI (2917)  
2900 GGTAATAGCGATGACTAATACGTAGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGCATAATGCCAGGCGGGCCATTTACCGTCAATTG

NdeI (3022)  
3000 ACGTCAATAGGGGCGTACTTGGCATATGATACACTTGTACTGCCAAGTGGGCGTGTACCGTAAATACTCCACCCATTGACGTCAATGAAAGTCC  
3100 CTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGCTGTTGGCGGTACGACGCGGGCCATTTACCGTAAGTTATGTAAC

PacI (3208)  
PstI (3201)  
**SdaI (3200)**      **BspLU11I (3218)**

3200 GCCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCC  
3300 CTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTGCG  
3400 CTCTCTGTTCGACCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCAGCTGTAGGTATCTCAGT

ApaLI (3532)

3500 TCGGTGTAGGTCGTTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACC  
3600 CGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGCC  
3700 TAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAA  
3800 ACCACCGCTGGTAGCGGTGTTTTTTTTGTTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTACGGGGTCTG

EagI (3968)  
PacI (3948)    SwaI (3957)    **NotI (3967)**

3900 ACGCTCAGTGGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTA  
4000 CATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCCCAGTG  
4100 CAAGTGCAGGTGCCAGAACATTTCTCTATCGAA