

pFUSE-hIgG1e13-Fc1

Plasmid containing a human engineered IgG1 Fc region

Catalog # pfc1-hg1e13

For research use only

Version 20K04-MM

PRODUCT INFORMATION

Content:

- 20 µg of pFUSE-hIgG1e13-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 activity¹. 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

- **hIgG1e13-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 decreased FcRn binding affinity at pH 6.0 and enhanced antibody clearance^{2,3}. The engineered pFUSE-hIgG1e13-Fc1 contains a single amino acid substitution: I253A.
- **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1α (EF-1α) core promoter⁴ and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat⁵. 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 mRNA⁶.
- **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 transcription⁷.

1. Carter PJ, 2006. Potent antibody therapeutics by design. *Nat Rev Immunol*. 6, 343-357.
2. Petkova SB, et al. (2006). Enhanced half-life of genetically engineered human IgG1 antibodies in a humanized FcRn mouse model: potential application in humorally mediated autoimmune disease. *Int Immunol* 18(12):1759-69. 3. Qiao S-W, et al., 2008. Dependence of antibody-mediated presentation of antigen on FcRnPNAS 105(27): 9337-9342. 4. 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. 5. 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. 6. 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. 7. 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](#)

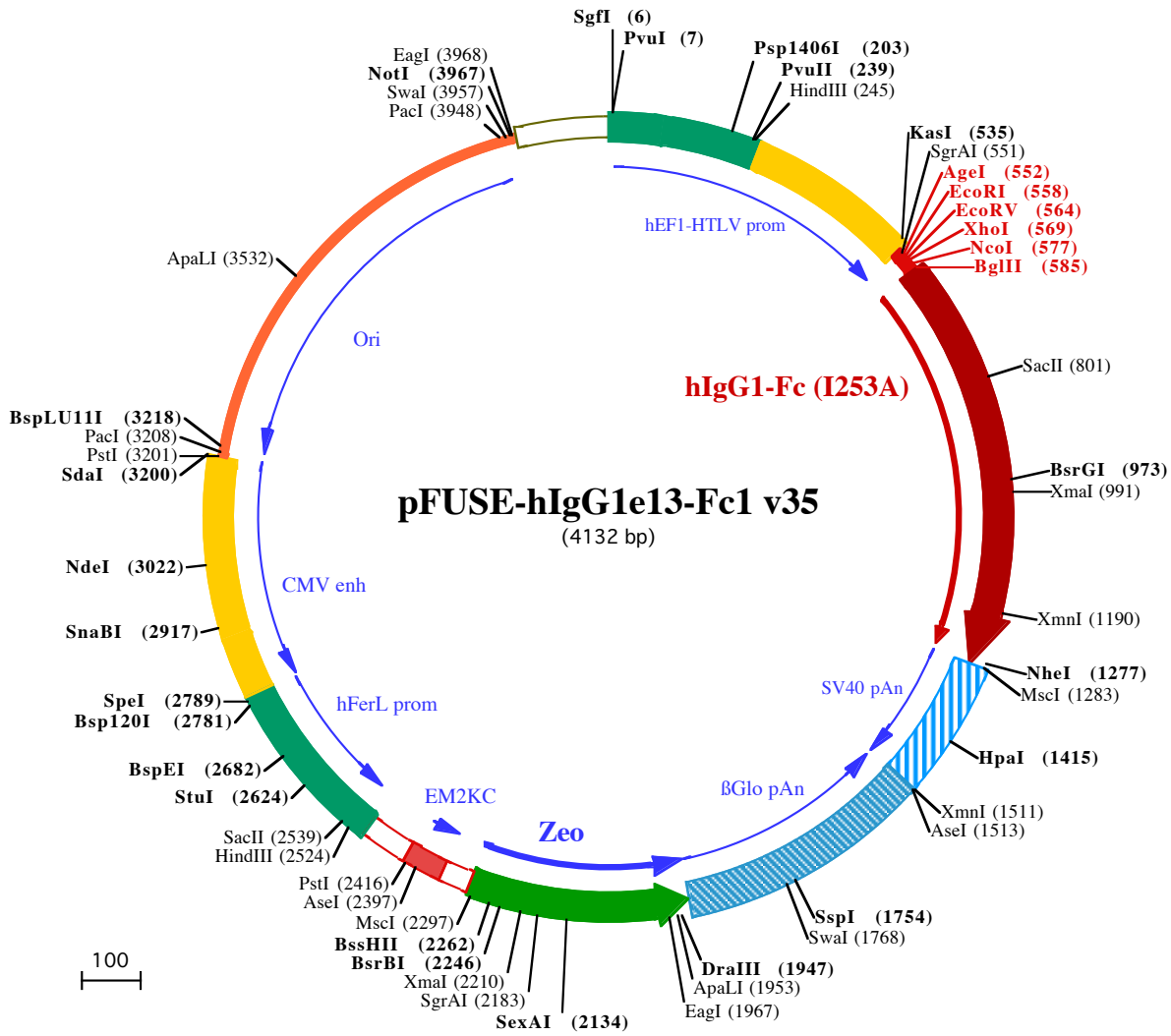
InvivoGen USA (Toll-Free): 888-457-5873

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



PvuI (7)
SgfI (6)
 1 GGATCTGCGATCGCTCCGGTCCCGTCAGTGGGACAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGTGGCAATTGAACGGTGCCTA
 101 GAGAAAGTGGCGGGGTAAGTGGAAAGTGATGTCGTGACTGGCTCCGCCCTTTTCCCGAGGGTGGGGGAGAACGTATATAAGTGCAGTAGTCGCC

HindIII (245)
Psp1406I (203) **PvuII (239)**
 201 GTGAACGTTCTTTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTCGAGGGGTCGCATCTCTCTTCACGCCGCCGCCCTACCTGAGGCC
 301 GCCATCCACGCCGGTTGAGTCGCTTCTGCCGCTCCCGCTGTGGTGCCTCTGAAGTGGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGTCCGAGACC
 401 GGGCCTTTGTCCGGCGCTCCCTTGAGGCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCCTGACCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

EcoRI (558) **KasI (535)** **AgeI (552)** **XhoI (569)** **BglII (585)**
SgrAI (551) **EcoRV (564)** **NcoI (577)**
 501 TCTGTTCTGCCCGTTACAGATCCAAGCTGTGACCGGCCCTACCTGAGATCACCGGTAATTCGATATCTCGAGCACCATGGTTAGATCTGACAAAAC
 601 CACACATGCCACCCTGCCAGCACCTGAAGTCTGGGGGACCGTCACTTCTCTTCCCCAAAACCAAGGACACCTCATGGCTCCCGGACCC
 4 H T C P P C P A P E L L G G P S V F L F P P K P K D T L M A S R T
 701 CTGAGGTCACATGCTGGTGGTGGACGTGAGCCACGAAGACCTGAGGTCAAGTCAACTGGTACGTGGACGGCGTGGAGGTGCATAATGCCAAGACAAA
 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 K
SacII (801)
 801 GCCGGGAGGAGCAGTACAACAGCACGTACCCTGTGGTCAAGCTCCTACCCTGCTGACCCAGGACTGGCTGAATGGCAAGGAGTACAAGTGAAGGTC
 70 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) **XmaI (991)**
 901 TCCAACAAGCCCTCCAGCCCCATCGAGAAAACCATCTCAAAGCCAAAGGGCAGCCCCGAGAACACAGGTGTACACCTGCCCCATCCGGGAGG
 104 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 T
 1001 AGATGACCAAGAACAGGTGACCTGACCTGCCTGGTCAAAGCTTCTATCCGAGCACATCGCCGTGGAGTGGGAGAGCAATGGCGCCGAGAACAA
 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 N
XmnI (1190)
 1101 CTACAAGACCAGCTCCCGTGGTGGTCCGACGGCTCTTCTCTCTACAGCAAGCTCACCGTGGACAAGAGCAGGTGGCAGCAGGGAAAGCTTCTC
 170 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)
 1201 TCATGCTCCGTGATGCACGAGGCTCTGCACAACCACTACACGCAGAAGGCTCTCCCTGTCTCCGGGTAATGAGTGTAGTGGCCAGACATGATAAG
 204 S C S V M H E A L H N H Y T Q K S L S L S P G K •
 1301 ATACATTGATGAGTTTGGACAAACCAACTAGAAATGCAGTGAATAAATGCTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATA

HpaI (1415)
 1401 AGCTGCAATAAACAAGTTAAACAACAATTCATTCTTTATGTTTCAGGTTCAAGGGGAGGTGGGGAGGTTTTTAAAGCAAGTAAAACCTCTACA

AseI (1513) **XmnI (1511)**
 1501 AATGTGGTATGGAATTAATTCTAAATACAGCATAGCAAAACTTAACTCCAATCAAGCCTCTACTTGAATCTTTTCTGAGGGATGAATAAGGCATA
 1601 GGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTCATGGAGTTTAAAGATATAGTGATTTTTCCCAAGGTTTGAACATAGCT

SspI (1754) **Swal (1768)**
 1701 CTTCAATTTCTTTATGTTTTAAATGCACGTGACCTCCACATTCCTTTTATAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATG
 1801 TTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTGTGGACTTAGGGAACAAAGAACCTTAAATAGAAATTGGA

ApaLI (1953) **DraIII (1947)** **EagI (1967)**
 1901 CAGCAAGAAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCTCTGCCCACAAAGTGCACGCAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCCCGCCCC
 125 D Q E E A V F H V C N G A P D R L A F E R G W
 2001 ACGGCTGCTCGCGATCTCGGTGATGCGCGGCCGAGGCGTCCCGAAGTTCGTGGACACGACCTCCGACCACTCGCGTACAGCTCGTCCAGGCCGG
 101 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)**
 2101 CACCACACCCAGGCCAGGGTGTGTCCGGCACCACTGGTCTGACCGCGCTGATGAACAGGGTCAAGTCCGCGTCCCGGACCAACCGCGAAGTCTGCC
 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
XmaI (2210) **BsrBI (2246)** **BssHII (2262)** **MscI (2297)**
 2201 TCCACGAAGTCCCGGGAGAACCAGCGGCTCGGTCCAGAAGTCCCGGCGACGTCGCGCGGTTGAGCACCAGGACCGGACTGGTCAACTTGG
 34 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 A
AseI (2397)
 2301 CCATGATGGCTCCTCctgtcaggagaggaagagaagaaggttagtacaattgCTATAGTGAAGTGTATTATACTATGCAGATATACTATGCCAATGATT
 14 M
PstI (2416)
 2401 AATTGTCAAAGTGGGCTGCAGggttcatagtgcacttttctgcactgccccatctcctgccaccctttccagcatagacagtcagtgacttacC

HindIII (2524) **SacII (2539)**
 2501 AAACCTCACAGGAGGAGAAGGCAGAAGCTTGAGACAGACCCGCGGACCGCCGAAGTGCAGAGGGACGTGGCTAGGGCGGCTTCTTTATGTTGCCCGG

StuI (2624) **BspEI (2682)**
 2601 CCCTCGAGGAGGGCGCTCGGGAGCCATGCGGCTGCGAGGAGGGGGCCGAGGCGGCGTGCCTGACCAATCCGGAGCAGATAGGAGT

SpeI (2789) **Bsp120I (2781)**
 2701 CTCAGCCCCCGCCCAAGCAAGGGGAAGTACGCGCTGTAGCGCCAGCGTGTGTGAAATGGGGCTTGGGGGGTGGGGCCCTGACTAGTCAAAA
 2801 CAAACTCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCGTGAAGTCAAACCGCTATCCACGCCATTGATGACTGCCAAAACCGCATCATCATG

SnaBI (2917)
 2901 GTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCCATAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTTACCGTCAATTGA

NdeI (3022)
 3001 CGTCAATAGGGGGCTACTTGGCATATGATACACTTGTACTGCAAGTGGGCGAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCC
 3101 TATTGGCGTACTATGGGAACATACGTCAATTATTGACGTCAATGGCGGGGGTCTTGGGGGTGAGCCAGGCGGGCCATTTACCGTAAAGTTATGTAACG

Pacl (3208)
PstI (3201)
SdaI (3200) BspLU11I (3218)
3201 CCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCTTGCTGGCGTTTTCCATAGGCTCCGCCCC
3301 TGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGCGTTTCCCCCTGGAAGCTCCCTCGTGCC
3401 TCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCCTTCTCCCTTCGGGAAGCGTGGCGCTTTCATAGCTCACGCTGTAGGTATCTCAGTT
ApaLI (3532)
3501 CGGTGTAGGTCGTTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTTCAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCC
3601 GGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCT
3701 AACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAA
3801 CCACCGCTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTACGGGGTCTGA
EagI (3968)
Pacl (3948) SwaI (3957) NotI (3967)
3901 CGCTCAGTGGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTAC
4001 ATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCCCAGTGC
4101 AAGTGCAGGTGCCAGAACATTTCTCTATCGAA