

pFUSE-hIgG1e12-Fc1

Plasmid containing a human engineered IgG1 Fc region

Catalog # pfc1-hg1e12

For research use only

Version 20K04-MM

PRODUCT INFORMATION

Content:

- 20 µg of pFUSE-hIgG1e12-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

- **hIgG1e12-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 6.0². The engineered pFUSE-hIgG1e12-Fc1 contains two amino acid substitutions: M428L and N434A.
- **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1α (EF-1α) core promoter³ and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat⁴. 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 mRNA⁵.
- **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 transcription⁶.

1. Carter PJ., 2006. Potent antibody therapeutics by design. Nature Reviews Immunology. Nat Rev Immunol. 6, 343-357.
2. Yeung Y.A. *et al.*, 2009. Engineering Human IgG1 Affinity to Human Neonatal Fc Receptor: Impact of Affinity Improvement on Pharmacokinetics in Primates. J. Immunol. 182: 7663-7671.
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



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

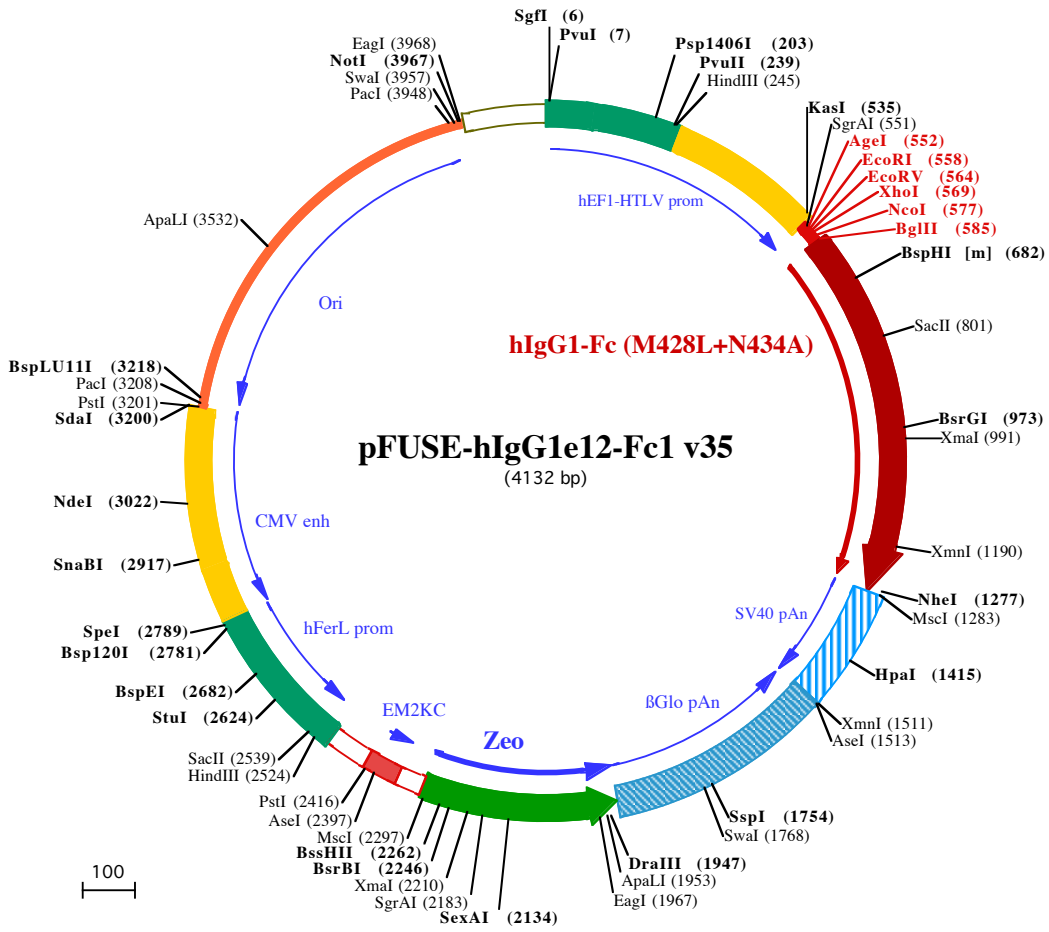
InvivoGen USA (Toll-Free): 888-457-5873

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

HindIII (245)
201 GTGAACGTTCTTTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTCGAGGGGTCGCATCTCTCTTACGCGCCCGCCCTACCTGAGGCC

Psp1406I (203) PvuII (239)
301 GCCATCCACGCCGGTTGAGTCGCTTCTGCCGCTCCCGCTGTGGTGCCTCTGAAGTGGTCCGCGCTAGGTAAGTTTAAAGCTCAGTGCAGAGCC

401 GGGCCTTTGTCCGGCGCTCCCTTGAGGCTACCTAGACTCAGCCGGCTCTCCACGCTTGCCTGACCTGCTTGTCAACTCTAGCTCTTTGTTTCGTTT

EcoRI (558) AgeI (552) XhoI (569) BglII (585)
KasI (535) SgrAI (551) EcoRV (564) NcoI (577)
501 TCTGTTCTGCCCGTTACAGATCCAAGCTGTGACCGGCCCTACCTGAGATCACCGGTAATTCGATATCTCGACACCATTGGTTAGATCTGACAAAAC

BspHI [m] (682)
601 CACACATGCCACCGTCCCGACCTGAAGTCTGGGGGACCGTCACTTCTCTTCCCCAAAACCAAGGACACCCTCATGATCTCCCGGAGCC

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 I S R T
701 CTGAGGTCACATCGTGGTGGTGGACGTGAGCCACGAGACCTGAGGTCAAGTCAACTGGTACGTGGACGGCTGGAGGTGCATAATGCCAAGACAAA

37 P E V T C 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
801 GCCGCGGAGGAGCAGTACAACAGCACGTACCGTGTGGTCAAGCTCCTCACCGTCTGCACAGGACTGGCTGAATGGCAAGGAGTACAAGTCAAGGTC

SacII (801)
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 TCCAACAAGCCCTCCGACGCCCCATCGAAGAAACATCTCAAAGCCAAAGGGCAGCCCCGAGAACCACAGGTGTACACCCTGCCCCATCCCGGGAGG

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 E
1001 AGATGACCAAGAACCAGGTGAGCTGACCTGCCTGGTCAAAGGCTTCTATCCAGCAGCATCGCCGTGGAGTGGGAGAGCAATGGGAGCCGGAGAACAA

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
1101 CTACAAGACCAGCTCCCGTGGACTCCGACGGCTCTTCTCTCTACAGCAAGCTCACCGTGGACAAGAGCAGGTGGCAGCAGGGGAACGTTCTTC

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
NheI (1277)
1201 TCATGCTCCGTGCTGCATGAGGCTCTGCACGCCACTACAGCAGAAGAGCCTCTCCCTGTCTCCGGTAAATGAGTGTAGCTGCTGCGCAGACATGATAAG

204 S C S V L H E A L H A H Y T Q K S L S P G K •
1301 ATACATTGATGAGTTGGACAACCAACTAGAATGCAAGTGAAGAAATGCTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATA

HpaI (1415)
1401 AGCTGCAATAAACAAGTTAAACAACAATTCGATTCTTTTATGTTTCAGGTTCCAGGGGAGGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACA

AseI (1513) XmnI (1511)
1501 AATGTGGTATGGAATTAATCTAAATACAGCATGAAAACTTAACTCCAAATCAAGCCTCTACTGAAATCCTTTCTGAGGATGAATAAGGCATA

1601 GGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTGACGCTCACCTTCTTCATGGAGTTTAAAGATATAGTGTATTTTCCCAAGTTTTGAACCTAGCT

SspI (1754) SmaI (1768)
1701 CTTCAATTTCTTATGTTTTAAATGCATGACCTCCACATTCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATG

1801 TTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGA

ApaLI (1953) DraIII (1947) EagI (1967)
1901 CAGCAAGAAAGCGAGCTTCTAGTCTTCTCAGTCTGCTCCTCTGCCACAAAGTGCACGAGTTGCCGGCCGGTGCAGGCGAACTCCCGCCCC

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 ACGGCTGCTCGCCGATCTCGGTATGCGCCGCGGAGGCGTCCCGAAGTTCGTGGACACGACCTCCGACCTCGGCGTACAGCTCGTCCAGGCCGG

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
2101 CACCCACACCCAGGCCAGGTTGTGTCGGCACCACCTGCTGACCCGCGCTGATGAACAGGGTCAAGTCTGCTCCCGACACACCCGGCAAGTCTGTC

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
2201 TCCACGAAGTCCCGGAGAACCCGAGCGGTGCTCCAGAACTGACCGCTCCGCGCAGCTCGCGCGGTGAGCACCAGGACCGTGGTCAACTGG

XmaI (2210) BsrBI (2246) BssHII (2262) MseI (2297)
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
2301 CCATGATGGCTCCTcgtcaggagaggaagagagagaggttagtacaattgTATAGTGGTGTATTATACTATGCAGATATACTATGCCAATGATT

1 M
2401 AATTGTCAAAGTGGGCTGCAgggttcatagtgcacttttctgcactgccccatctcctgccaccctttccaggcatagacagtcagtgacttacC

HindIII (2524) SacII (2539)
2501 AAATCAGAGGAGGAGAAGGAGAAGCTTGAGACAGACCCGCGGACCGCGAACTGCGAGGGGACGTGGCTAGGGCGGCTCTTTTATGGTGCAGCCG

StuI (2624) BspEI (2682)
2601 CCCTCGGAGGAGGCGCTCGGGAGGCTAGCGGCAATCTGCGGTGGCAGGAGCGGGCCGAAGCCGTGCCTGACCAATCCGGAGCACATAGGAGT

Bsp120I (2781) SpeI (2789)
2701 CTCAGCCCCCGCCCCAAAGCAAGGGGAAGTCAAGCGCTGTAGCGCCAGCGTGTGTGAAATGGGGCTTGGGGGGTGGGGCCCTGACTAGTCAAAA

2801 CAAACTCCATTGACGTCATGGGGTGGAGACTTGGAAATCCCGTGGTCAAACCGTATCCACGCCATTGATGACTGCCAAAACCGCATCATCATG

SnaBI (2917)
2901 GTAATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCATAAGGTCACTGACTGGGCATAATGCCAGCGGGCCATTACCGTCATTGA

NdeI (3022)
3001 CGTCAATAGGGGCGTACTTGGCATATGATACACTTGTACTGCCAAGTGGGCAAGTTTACCCTAAATCTCCACCCATTGACGTCATGGAAAGTCCC

3101 TATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGGCGGGGGTCGTTGGCGGTTCAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACG

PacI (3208)

PstI (3201)

SdaI (3200)

BspLU11I (3218)

3201 CCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGAACCGTAAAAAGCCGCGTTGCTGGCGTTTTCCATAGGCTCCGCCCC

3301 TGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCTGGAAGCTCCCTCGTGCC

3401 TCTCCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCCTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTT

ApaLI (3532)

3501 CGGTGTAGGTCGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTACGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCC

3601 GGTAAGACACGACTTATGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCT

3701 AACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAA

3801 CCACCGCTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTACGGGGTCTGA

EagI (3968)

PacI (3948) SwaI (3957)

NotI (3967)

3901 CGCTCAGTGGAAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTAC

4001 ATCTGTGTGTTGGTTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCCGAGTGC

4101 AAGTGCAGGTGCCAGAACATTTCTCTATCGAA