

pFUSE-hIgG4-Fc1

Plasmid designed for the construction of Fc-Fusion proteins

Catalog # pfuse-hg40fc1

For research use only

Version 20K05-MM

PRODUCT INFORMATION

Content:

- 20 µg of pFUSE-hIgG4-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 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 interactions¹.
- **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1α (EF-1α) core promoter² and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat³. 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 mRNA⁴.
- **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 transcription⁵.

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

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

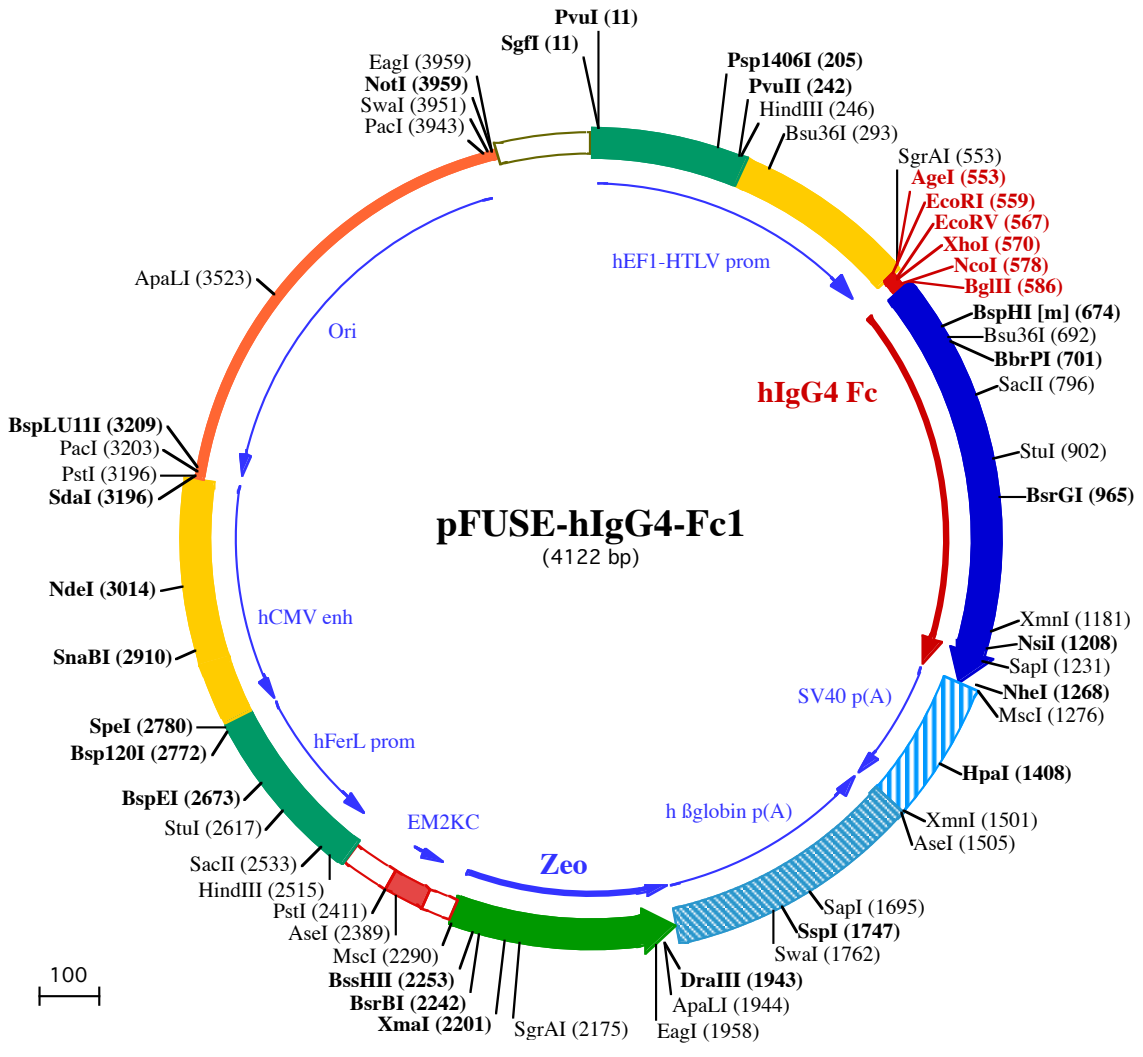
This antibiotic can be used for *E. coli* at 25 µg/ml in liquid or solid media and at 50-200 µg/ml to select Zeocin™-resistant mammalian cells.

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 (11)
SgfI (11)
 1 GGATCTGCATCGCTCCGGTGCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
 101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGTCTGTACTGGCTCCGCTTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (246)
Psp1406I (205) **PvuII (242)** **Bsu36I (293)**
 201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACCGCGCCGCCCTACCTGAGGGC
 301 GCCATCCACGCCGGTTGAGTCGCGTTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAACTGCGTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC
 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTCCTGACCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

EcoRI (559)
AgeI (553) **XhoI (570)**
SgrAI (553) **EcoRV (567)** **NcoI (578)** **BglII (586)**
 501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTGAATTCGATATCTCGAGCACCATTGATCTCCCCCATGC
 1 P P C
BbrPI (701)

BspHI [m] (674) **Bsu36I (692)**
 601 CCATCATGCCAGCACCTGAGTTCCTGGGGGACCATCAGTCTTCTGTTCCCCCAAACCAAGGACACTCTCATGATCTCCGGACCCCTGAGGTCA
 4 P S C P A P E F L G G P S V F L F P P K P K D T L M I S R T P E V
SacII (796)
 701 CGTGCCTGGTGGTGGACGTGAGCCAGGAAGACCCCGAGGTCCAGTCAACTGGTACGTGGATGGCGTGGAGGTGCATAATGCCAAGACAAAGCCGCGGGA
 37 T C V V V D V S Q E D P E V Q F N W Y V D G V E V H N A K T K P R E
StuI (902)
 801 GGAGCAGTTCAACAGCACGTACCGTGTGGTCAGCGTCTCACCCTGCACCAGGACTGGCTGAACGGCAAGGAGTACAAGTGAAGGTCTCCAACAAA
 70 E Q F 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 S N K
BsrGI (965)
 901 GGCTCCCGTCTCCATCGAGAAAACCTCTCAAAGCAAAGGGCAGCCCCGAGAGCCACAGGTGTACACCCTGCCCCATCCAGGAGGAGATGACCA
 104 G L P S S I E K T I S K A K G Q P R E P Q V Y T L P P S Q E E M T
 1001 AGAACCAGTCACTGACCTGCTGGTCAAAGCTTCTACCCAGCGACATCGCCGTGGAGTGGGAGAGCAATGGGAGCCGAGACAACCTACAAGAC
 137 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 Y K T
XmnI (1181)
 1101 CACGCCTCCCGTGTGGACTCCGACGGCTCCTTCTCTCTACAGCAGCTAACCGTGGACAAGAGCAGGTGGCAGGAGGGAATGTCTTCTCATGTCC
 170 T P P V L D S D G S F F L Y S R L T V D K S R W Q E G N V F S C S

NsiI (1208) **SapI (1231)** **NheI (1268)** **MscI (1276)**
 1201 GTGATGCATGAGGCTCTGCACAACCACTACACACAGAAGACCTCTCCCTGTCTCTGGGTAATAAAGCTAGCTGGCCAGACATGATAAGATACATTGA
 204 V M H E A L H N H Y T Q K S L S L S L G K •
 1300 TGAGTTTGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAAT

HpaI (1408)
 1400 AAACAAGTTAAACAACAACAAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTA

AseI (1505) **XmnI (1501)**
 1500 TGAATTAATTCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGG
 1600 GGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTCATGGAGTTTAAAGATATAGTGTATTTTCCCAAGGTTTGAACCTAGCTCTTCATTTT

SspI (1747) **SwaI (1762)**
 1700 TTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTTAGTAAAATATTAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATT
 1800 AGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAAGGAACCTTAAATAGAAATTGGACAGCAAGAA

ApaLI (1944) **DraIII (1943)** **EagI (1958)**
 1900 AGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTGCCACAAAGTGCACGCAGTTGCCGGCCGGGTGCGCAGGGCGAACTCCCCCCCCACGGCTGCT
 125 D Q E E A V F H V C N G A P D R L A F E R G W P Q E
 2000 CGCCGATCTCGGTATGGCCGGCCGGAGGCGTCCCGAAGTTCGTGGACACGACCTCCGACCACTCGCGTACAGCTCGCCAGGCCGCGCACCCACAC
 98 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 V W V
SgrAI (2175)
 2100 CCAGGCCAGGGTGTGTCCGGCACCACTGGTCTGGACCGCGCTGATGAACAGGGTACGTCGTCGCCGACCAACCCGGCGAAGTGTCTCCACGAAG
 65 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 E V F

XmaI (2201) **BsrBI (2242)** **BssHII (2253)** **MscI (2290)**
 2200 TCCCGGGAGAACCCGAGCCGGTCCGCTCCAGAACTCGACCGCTCCGGCGACGTGCGCGCGGGTGGAGCACCGGAACGGCACTGGTCAACTTGGCCATGATGG
 31 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 M
AseI (2389)
 2300 CTCCTCctgtcaggagaggaagagaagaaggttagtacaattgCTATAGTGAGTTGTATTACTATGCAGATATACTATGCCAATGATTAATTGTCAA
PstI (2411)
 2400 ACTAGGGCTGCAGggttcatagtgccacttttctgcactgccccatctctgcccaccctttcccaggcatagacagtcagtgacttacCAAACCTACA
HindIII (2515) **SacII (2533)**
 2500 GGAGGGAGAAGGCAGAAGCTTGAGACAGACCCGCGGGACCGCCGAACCTGCGAGGGGACGTGGCTAGGGCGGCTTCTTTTATGGTGCGCCGGCCCTCGGAG
StuI (2617) **BspEI (2673)**
 2600 GCAGGGCGCTCGGGGAGCCTAGCGCCAATCTGCGGTGGCAGGAGCGGGGCCGAAGGCCGTGCCTGACCAATCCGGAGCACATAGGAGTCTCAGCCCC
SpeI (2780) **Bsp120I (2772)**
 2700 CCGCCCCAAGCAAGGGGAAGTCACGCGCTGTAGCGCCAGCGTGTGTGAAATGGGGGCTTGGGGGGTGGGGCCCTGACTAGTCAAAACAAACTCCC
 2800 ATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCG
SnaBI (2910)
 2900 ATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCCATAAGGTCATGTACTGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAG
NdeI (3014)
 3000 GGGGCGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGCAGTTTACCGTAAATACTCCACCATTGACGTCAATGGAAAGTCCCTATTGGCGT
PacI (3203) **PstI (3196)** **SdaI (3196)**
 3100 TACTATGGGAACATACGTCAATTATTGACGTCAATGGCGGGGGTCTTGGCGGTGAGCCAGGCGGGCCATTTACCGTAAAGTTATGTAACGCCCTGCAGGT
BspLU11I (3209)
 3200 TAATTAAGAACATGTGAGCAAAAGGCCAGAAAAGGCCAGAACCGTAAAAAGCCGCTTGTGGCCTTTTTCCATAGGCTCCGCCCCCTGACGAGCA
 3300 TCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTT
 3400 CCGACCCTGCCGTTACCGGATACCTGTCCGCCTTCTCCCTTCGGGAAGCGTGGCGCTTTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGG
ApaLI (3523)
 3500 TCGTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGTAGTCCAACCCGGTAAGACA
 3600 CGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGC
 3700 TACTACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTG
 3800 GTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTG
EagI (3959) **PacI (3943)** **Swal (3951)** **NotI (3959)**
 3900 GAACGAAAACCTACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAAATATCTTTATTTTATTACATCTGTGTG
 4000 TTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAAACTAGCAAATAGGCTGTCCCCAGTGCAAGTGCAGG
 4100 TGCCAGAACATTTCTCTATCGAA