

pFUSE-rtIgG2B-Fc1

Plasmid designed for the construction of rat IgG2B Fc-Fusion proteins

Catalog # pfuse-rtg2bfc1

For research use only

Version 20K05-MM

PRODUCT INFORMATION

Content:

- 20 µg of pFUSE-rtIgG2B-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 a sequence encoding a given protein to the Fc region of an immunoglobulin.

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. These cells 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.

PLASMID FEATURES

- **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1α (EF-1α) core promoter¹ and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat². 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.
5' - Age I, Eco RI, Eco RV, Xho I, Nco I, Bgl II - 3'
- **rtIgG2B Fc (rat):** The Fc region comprises the CH2 and CH3 domains of the IgG2B 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.

- **SV40 pAn:** the Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA³.
- **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 transcription⁴.

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.

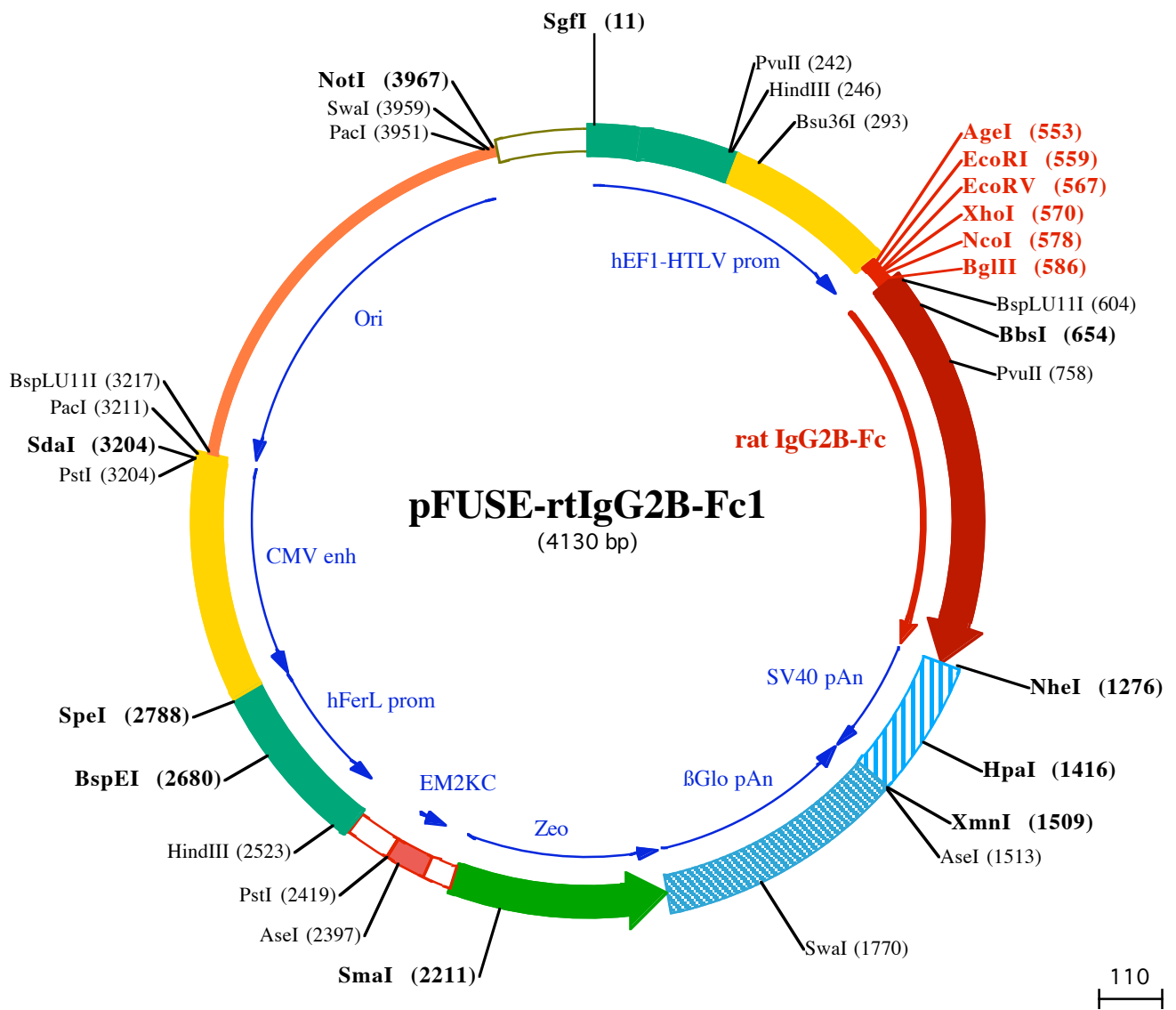
References:

1. Kim, D.W. *et al.* (1990). *Gene* 2: 217-223.
2. Takebe, Y. *et al.* (1988). *Mol. Cell Biol.* 1: 466-472.
3. Carswell, S., and Alwine, J.C. (1989). *Mol. Cell Biol.* 10: 4248-4258.
4. Yu J & Russell JE. (2001). *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





SgfI (11)
1 GGATCTGCGATCGCTCCGGTCCCGTCAGTGGCGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGTTCGCAATTGAACGGGTGCCTA
101 GAGAAGGTGGCGGGGTAACCTGGGAAAGTGATGTCGTGTACTGGCTCCGCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (246)
PvuII (242)
201 GTGAACGTTCTTTTTCGCAACGGGTTTGC CGCAGAAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACCGCGCCCGCCCTACCTGAGGCC
301 GCCATCCACGCGGGTTGAGTCGCGTTCTGCCCTCCCGCTGTGGTGCCTCTGAAGCTCGCTCCGCGCTAGGTAAGTTTAAAGCTCAGGTCGAGACC
401 GGGCTTTGTCCGGCTCCCTTGAGCTACCTAGACTCAGCGGCTCCACGCTTTCCTGACCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

EcoRI (559) XhoI (570) BglII (586)
AgeI (553) EcoRV (567) NcoI (578)
501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGGCGCTACCTGAGATCACCGTGAATTCGATATCTCGAGCACCATGGTTAGATCTCTACATGCT
1►ProThr Cys

BspLU111 (604)
601 CCTACATGTCAAAATGCCAGTTCCTGAACTCTGGTGGACCATCTGCTTCATCTCCCGCAAAGCCCAAGGACATCTCTTGTATCTCCAGAACG
4►ProThr CysHisLysCysProValProGluLeuLeuGlyGlyProSerValPheIlePheProProLysProLysAspIleLeuLeuIleSerGlnAsnA

BbsI (654)
PvuII (758)
701 CCAAGGTCACGTGTGGTGGTGGATGTGAGCGAGGAGCGGACGCTCCAGTTCAGCTGGTTTGTGAACAACGTAGAAGTACACAGCTCAGACACA
37►IaLysValThrCysValValValAspValSerGluGluGluProAspValGlnPheSerTrpPheValAsnAsnValGluValHisThrAlaGlnThrGln
801 ACCCGTGAGGAGCAGTACAACAGCACCTTACAGTGGTTCAGTGCCTCCCATCCAGCACGAGGACTGGATGAGCGCAAGGAGTTCAAATGCAAGGTC
70►nProArgGluGluGlnTyrAsnSerThrPheArgValValSerAlaLeuProIleGlnHisGlnAspTrpMetSerGlyLysGluPheLysCysLysVal
901 AACAAACAAAGCCCTCCAAAGCCCATCGAAGAACCATCTCAAACCCAAAGGCTAGTCAGAAACACAGGTATACGTCATGGTCCACCGCAGAGC
104►AsnAsnLysAlaLeuProSerProIleGluLysThrIleSerLysProLysGlyLeuValArgLysProGluValTyrValMetGlyProProThrGluG
1001 AGTTGACTGAGCAAACGGTTCAGTTGACCTGCTGACCTCAGGCTTCCTCCATACGACATCGGTGTGGAGTGAGCAGCAACGGGCATATAGAAAAGAA
137►InLeuThrGluGlnThrValSerLeuThrCysLeuThrSerGlyPheLeuProAsnAspIleGlyValGluTyrThrSerAsnGlyHisIleGluLysAs
1101 CTACAAGAACACCGAGCCAGTGTGACTCTGACGGTCTTTCTCATGTACAGCAAAGCTCAATGTGAAAGGAGCAGGTGGGATAGCAGAGCGCCCTTC
170►nTyrLysAsnThrGluProValMetAspSerAspGlySerPhePheMetTyrSerLysLeuAsnValGluArgSerArgTrpAspSerArgAlaProPhe

NheI (1276)
1201 GTCTGCTCCGTGTCACGAGGGTCTGCACAATCACCGTGGAGAAGAGCATCTCCCGGCTCCGGTAAATGAGCTAGCTGACAGCATGATAAGAT
204►ValCysSerValValHisGluGlyLeuHisAsnHisHisValGluLysSerIleSerArgProProGlyLys●●●

1301 ACATTGATGAGTTGGCAAACCAACTAGAATGCAGTGAATAAATGCTTTATTTGTGAATTTGTGATGCTATTGCTTTATTTGTAACCATATAAG

HpaI (1416)
1401 CTGCAATAAACAAGTTAAACAACAACAATTGCATTCATTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAA

AseI (1513)
XmnI (1509)
1501 TGTGGTATGGAAATTAATTTCAAATACAGCATAGCAAACTTTAACTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGG
1601 CATCAGGGGCTGTTGCAATGTGCATTAGCTGTTTGCAGCCTCACCTCTTTCATGGAGTTAAGATATAGTGATTTTTCCCAAGGTTTGAAGTCTCT

SwaI (1770)
1701 TCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTATGATAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTT
1801 TTTTATTAGGCAGAAATCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAAGGAACCTTAAATAGAAATGGACA

1901 GCAAGAAAGCGAGCTTCTAGCTTATCTCAGTCTGCTCCTTGCACAAAGTGCACGCAGTTGCGGCCGGTTCGCGCAGGGCGAAGTCCCGCCCCAC
125►●●●AspGlnGluGluAlaValPheHisValCysAsnGlyAlaProAspArgLeuAlaPheGluArgGlyTrpP
2001 GGCTGCTCGCGATCTCGGTCATGGCCGGCCGGAGGCTCCCGAAGTCTGTTGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCCCGCA
100►roGlnGluGlyIleGluThrMetAlaProGlySerAlaAspArgPheAsnThrSerValValGluSerTrpGluAlaTyrLeuGluAspLeuGlyArgVa
2101 CCCACACCCAGGCCAGGGTGTGTCGGCACCACTGTTCTGACCGCGTGAACAGGGTACGTCGTCGGGACACACCGGCAAGTCTGCTCCTC
67►I TrpValTrpAlaLeuThrAsnAspProValValGlnAspGlnValAlaSerIlePheLeuThrValAspAspArgValValGlyAlaPheAspAspGlu

SmaI (2211)
2201 CACGAAGTCCCGGGAGAACCCGAGCCGGTCCGAGAACTCGACCGCTCCGGCGAGCTGCGCGCGGTGAGCACCGGAACGGCACTGGTCAACTTGGCC
34►ValPheAspArgSerPheGlyLeuArgAspThrTrpPheGluValAlaGlyAlaValAspArgAlaThrLeuValProValAlaSerThrLeuLysAlaM
2301 ATGATGGCTCCTCctgtcaggagaggaaagagaagaaggttagtacaattgCTATAGTGAGTTGATTATACTATGAGATATACTATGCCAATGATTAA
0►et

PstI (2419)
2401 TTGTCAAACAGGGCTGCAgggttcatagtgccacttttctgactgccccatctctgccccacctttccaggcatagacagtcagtgacttacCAA

HindIII (2523)
2501 ACTCACAGGAGGGAGAAGGCTTGAGACAGACCCCGGGACCGCGAACTGCGAGGGGAGCTGGCTAGGGCGGCTCTTTTATGGTGC CGCGGCC

BspEI (2680)
2601 CTCGGAGGCAGGGCTCGGGGAGGCTAGCGGCAATCTGCGGTGGCAGGAGGGGGCCGAAGGCCGTGCCTGACCAATCCGGAGCACATAGGAGTCT

SpeI (2788)
2701 CAGCCCCCGCCCAAAGCAAGGGGAAGTACGCGCTGTAGCGCCAGCGTGTGTGAAATGGGGCTTGGGGGGTGGGGCCCTGACTAGTCAAAAACA
2801 AACTCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGT
2901 AATAGCGATGACTAATACGTAGTGTACTGCCAAGTAGAAAGTCCATAAGGTCATGTACTGGCATAATGCCAGCGGGCCATTTACCGTCATTGACC
3001 TCAATAGGGGCGTACTTGGCATATGATACACTTGTACTGCAAGTGGCGAGTTTACCGTAAATACTCCACCCATTGACGTCAATGAAAAGTCCCTAA

PstI (3204)
SdaI (3204)
3101 TTGGCGTTACTATGGAAACATACGTCATTATTGACGTCAATGGCGGGGCTGTTGGCGGTACAGCAGCGGGCCATTTACCGTAAGTTATGTAACGCC

PacI (3211) BspLU11I (3217)

3201 TGCAGGTTAATTAAAGACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTCCATAGGCTCCGCCCCCTG
3301 ACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCCTGGAAGCTCCCTCGTGCCTC
3401 TCCTGTTCCGACCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTG
3501 GTGTAGGTCGTTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGG
3601 TAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCTAA
3701 CTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACC
3801 ACCGCTGGTAGCGGTGGTTTTTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACG

PacI (3951) SwaI (3959) NotI (3967)

3901 CTCACTGGAAACGAAAACCTCACGTTAAGGGATTTTGGTCAAGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAAATATCTTTATTTTCATTACAT
4001 CTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCCAGTGCAA
4101 GTGCAGGTGCCAGAACATTTCTCTATCGAA