



## 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.

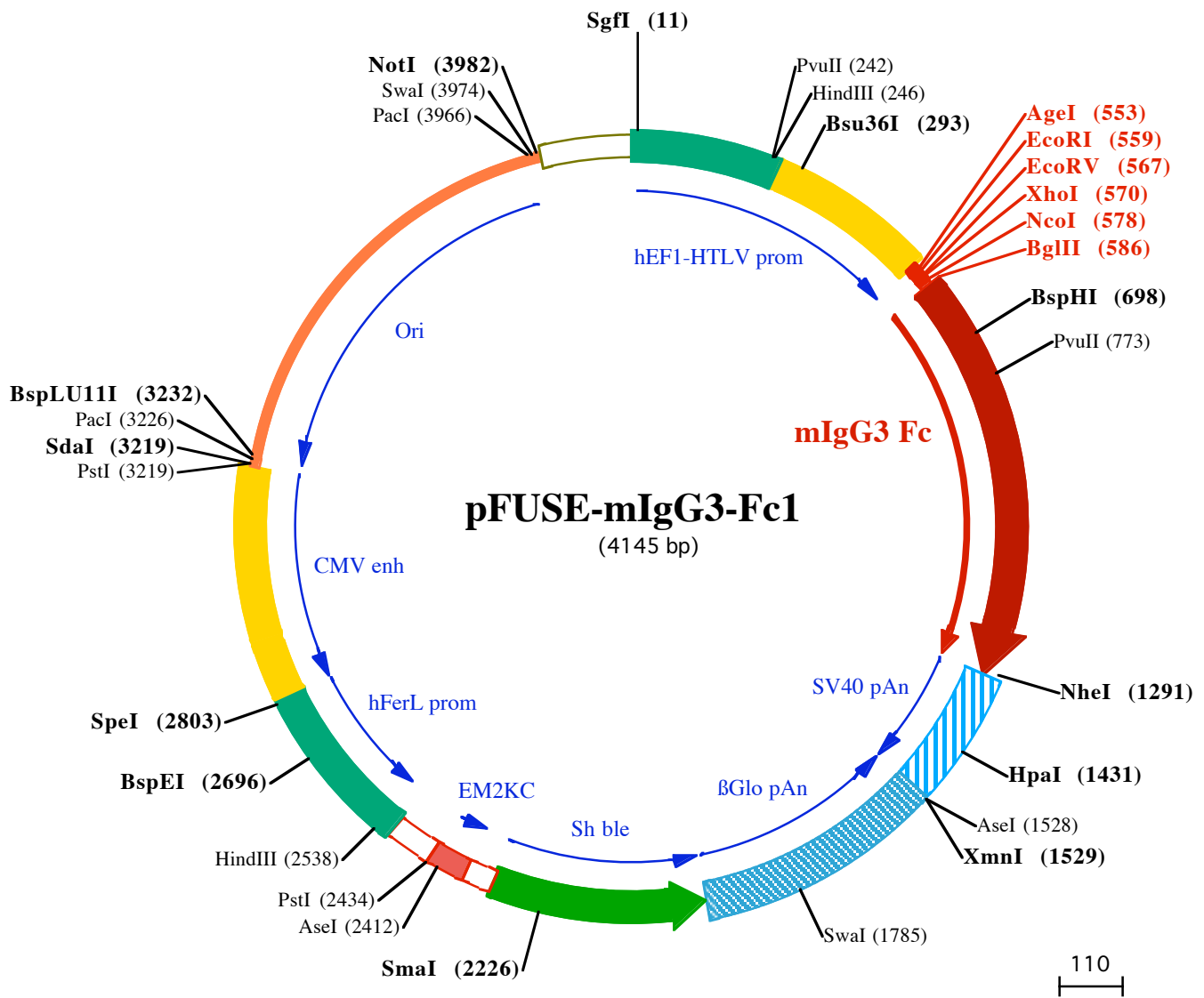
## RELATED PRODUCTS

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)



SgfI (11)  
1 GGATCTGCGATCGCTCCGGTGCCTGAGTGGGCGAGCGCACATCGCCACAGTCCCGAGAAGTTGGGGGAGGGTTCGCAATTGAACGGTGCCTA  
101 GAGAAGGTGGCGGGGTAACCTGGGAAAGTGATGTCGTGTACTGGCTCCGCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGGC

HindIII (246) PvuII (242) Bsu36I (293)  
201 GTGAACGTTCTTTTTGCAACGCGTTTGCCTCCGAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACGCGCCCGCCCTACCTGAGGCC  
301 GCCATCCACGCGGGTTGAGTGCCTTCTGCCCTCCCGCTGTGGTGCCTCTGAAGTCCGCTCCGCGCTAGGTAAGTTTAAAGCTCAGGTCGAGACC  
401 GGGCTTTGTCGGCGCTCCCTGGAGCTACCTAGACTCAGCGGCTCTCCACGCTTTCCTGACCTGCTTGTCTCAACTCTACGCTTTGTTTCGTTT

EcoRI (559) XhoI (570) BglII (586)  
AgeI (553) EcoRV (567) NcoI (578)  
501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCTACCTGAGATCACCGTGAATTCGATATCTCGAGCACCATGGTTAGTCTCTCTAGAATA  
1ProArgIle  
BspHI (698)  
601 CCCAAGCCAGTACCCCGAGTTCTTATGCCACCTGGTAACATCTTGGGTGGACCATCCGCTCTTATCTTCCCCCAAGCCCAAGGATGCACCTCA  
4ProLysProSerThrProProGlySerSerCysProProGlyAsnIleLeuGlyGlyProSerValPheIlePheProLysProLysAspAlaLeuM  
PvuII (773)  
701 TGATCTCCCTAACCCCAAGTTACGTGTGGTGGTGGATGTGAGCGAGGATGACCCAGATGTCCATGTGAGCTGGTTGTGGCAACAAAGAAGTACA  
377e t l eSerLeuThrProLysValThrCysValValValAspValSerGluAspAspProAspValHisValSerTrpPheValAspAsnLysGluValHis  
801 CACAGCTGGACACAGCCCGTGAAGCTCAGTACAACAGTACTTCCGAGTGGTTCAGTGCCTCCCATCCAGCACAGGACTGGATGAGGGCAAGGAG  
707sThrAlaTrpThrGlnProArgGluAlaGlnTyrAsnSerThrPheArgValValSerAlaLeuProIleGlnHisGlnAspTrpMetArgGlyLysGlu  
901 TTCAAATGCAAGGTCAACAACAAGCCCTCCAGCCCATCGAGAGAACCATCTCAAAACCCAAAGGAGAGCCAGACCTCAAGTATACACCATA  
104PheLysCysLysValAsnAsnLysAlaLeuProAlaProIleGluArgThrIleSerLysProLysGlyArgAlaGlnThrProGlnTyrThrIleP  
1001 CCCCACCTCGTGAACAATGTCCAAGAAGAAGTTAGTCTGACCTGCCTGGTCAACCACTTCTTCTGAAAGCCATCAGTGTGGAGTGGGAAAGAACGG  
137ProProProArgGluGlnMetSerLysLysLysValSerLeuThrCysLeuValThrAsnPhePheSerGluAlaIleSerValGluTrpGluArgAsnG  
1101 AGAAGTGGAGCAGGATTACAAGAACAACCTCCACCTCCTGGACTCAGATGGGACCTACTTCTCTACAGCAAGCTCACTGGGTGATACAGACAGTTGGTTG  
170yGluLeuGluGlnAspTyrLysAsnThrProProIleLeuAspSerAspGlyThrTyrPheLeuTyrSerLysLeuThrValAspThrAspSerTrpLeu  
NheI (1291)  
1201 CAAGGAGAAATTTTACCTGCTCCGTGGTGCATGAGGCTCTCCATAACCACACACAGAAGAACCCTGTCTCGCTCCCTGGTAAATGAGCTAGCTGGC  
204GlnGlyGluIlePheThrCysSerValValHisGluAlaLeuHisAsnHisHisThrGlnLysAsnLeuSerArgSerProGlyLys  
1301 CAGACATGATAAGATACATTGATGAGTTGGACAAACCACAACCTAGAATGCAGTGAAGAAATGCTTTATTTGTAAATTTGTGATGCTATTGCTTTATT

HpaI (1431)  
1401 TGTAACCATTATAAGCTGCAATAAACAAGTTAACACAACAATTGCATTTCATTTTATGTTTCAGGTTTCAGGGGAGGTTGGGGAGGTTTTTAAAGCAAG

AseI (1528) XmnI (1529)  
1501 TAAAACCTCTACAATGTGGTATGGAATTAATTCTAAAATACAGCATAGCAAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGA  
1601 TGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCCAAG

SwaI (1785)  
1701 GTTTGAACTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCA  
1801 ATGAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGCCCTTCAATAATCCCCAGTTTAGTAGTTGGACTTAGGAAACAAAGAACCTTT  
1901 AATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTATCTCAGTCTGCTCTGCCACAAGTGCACGAGTTGCGCGCCGGTTCGCGCAGGGCG  
125AspGlnGluGluAlaValPheHisValCysAsnGlyAlaProAspArgLeuAlaP  
2001 AACTCCCGCCCGCCGCTGCTCGCCGATCTCGGTTCATGGCCGCGCCGAGGCGTCCCGGAAGTTCTGGACACGACTCCGACCTCGCCGCTACAGCT  
105heGluArgGlyTrpProGlnGluGlyIleGluThrMetAlaProGlySerAlaAspArgPheAsnThrSerValValGluSerTrpGluAlaTyrLeuGlu  
2101 CGTCCAGGCGCCGACCCACACCCAGGCCAGGGTGTGTCGGCACCCTGCTGACCGCGCTGATGAACAGGGTACGTCGTCGGGACCCACCC  
72uAspLeuGlyArgValTrpValTrpAlaLeuThrAsnAspProValValGlnAspGlnValAlaSerIlePheLeuThrValAspAspArgValValGly  
SmaI (2226)  
2201 GCGAAGTCTCTCCACGAAGTCCCGGAGAACCCGAGCCGCTCGTCCAGAAGTCCGACCGCTCCGGCGACGTCGCGCGGGTGAACCCGGAACCGCA  
39AlaPheAspAspGluValPheAspArgSerPheGlyLeuArgAspThrTrpPheGluValAlaGlyAlaValAspArgAlaThrLeuValProValAlaSer  
2301 CTGGTCAACTTGCCATGATGGCTCCTCctgtcaggagaggaagaagaaggtagtacaattgCTATAGTGAGTTGATTATACTATGCAGATATAC  
5erThrLeuLysAlaMet  
AseI (2412) PstI (2434)  
2401 TATGCCAATGATTAATGTCAAACCTAGGGCTGCAgggttcatagtgccacttttctgactgccccatctctgccccctttccaggcatagacag  
HindIII (2538)  
2501 tcagtacttacAAAACCTCACAGGGGAGAAGCGAGAAGCTTGAGACAGCCCGGGACCGCGAACTGCGAGGGGACGTGGCTAGGGCGGCTCTTTT

BspEI (2696)  
2601 TATGGTCCCGCCCTCGGAGGCAGGGCGCTCGGGAGCCCTAGCGCCAACTGCGGTGGCAGGAGCGGGCCGAAGCCGCTGCTGACCAATCCGG  
2701 AGCACATAGGAGTCTCAGCCCCCGCCCAAGCAAGGGGAAGTACGCGCTGTAGCGCCAGCGTGTGTGAAATGGGGCTTGGGGGGTGGGGCCC

SpeI (2803)  
2801 TGACTAGTCAAAACAACTCCCATTCAGCTCAATGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCACGCCATTGATGACTGCCAAAA  
2901 CCGCATCATATGGTAATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCCATAAAGTTCATGACTGGGCATAATGCCAGGCGGGCCAT  
3001 TTACCGTCATTGACGTCATAGGGGGGCTACTTGGCATATGATACACTTGTACTGCCAAGTGGGCAGTTTACCCTAAATCTCCACCCATTGACGCT  
3101 AATGGAAGTCCCTATTGGCGTACTATGGAAACATACGTCATTATTGACGTCATAGGGCGGGGCTGTTGGCGGTGACGAGCGGGCCATTTACCGT

PacI (3226)

PstI (3219)

**SdaI (3219)**                      **BspLU11I (3232)**

3201 AAGTTATGTAACGCCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGTTGCTGGCGTTTTCCATA

3301 GGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAGATACCAGGCGTTTTCCCCCTGGAAG

3401 CTCCCTCGTGCCTCTCCTGTTCCGACCCCTGCCGCTTACCGGATACCTGTCGCGCTTTCTCCCTTCGGGAAGCGTGCGCTTTCTCATAGCTCACGCTGT

3501 AGGTATCTCAGTTCGGTGTAGGTCGTTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTTCAGCCCAGCCGCTGCGCTTATCCGGTAACTATCGTC

3601 TTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTT

3701 GAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGA

3801 TCCGCAAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTT

---

PacI (3966)    SmaI (3974)    **NotI (3982)**

3901 CTACGGGCTGACGCTCAGTGAACGAAACTCACGTTAAGGATTTTGGTCATGGCTAGTTAATTAACATTTAATCAGCGGCCCAATAAAATATCT

4001 TTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAATAGG

4101 CTGTCCCAGTGAAGTGCAGGTGCCAGAACATTTCTCTATCGAA