

METHODS

□□DDVPTG□□EIV□VSH□VDR□

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

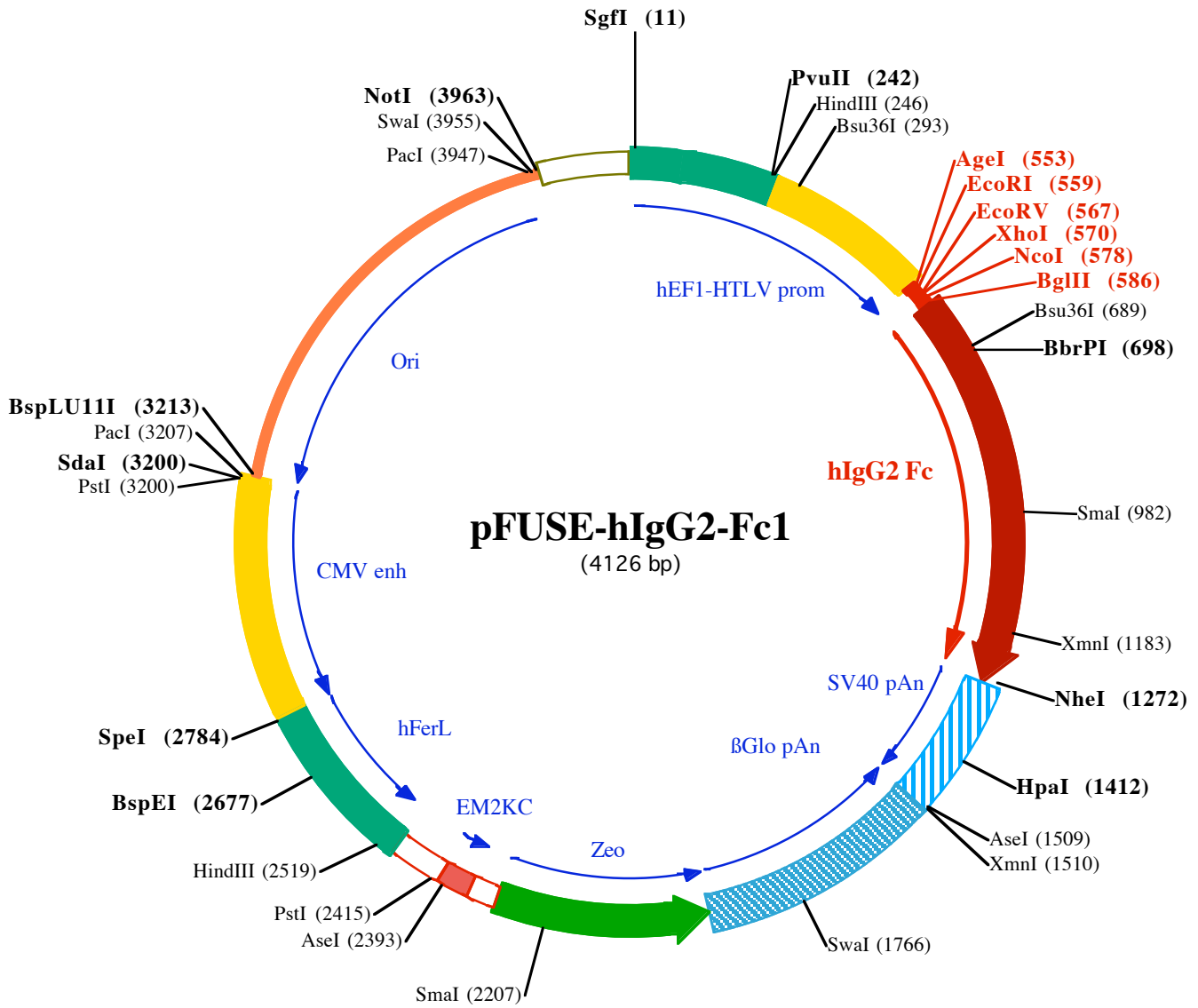
This antibiotic can be used for *E. coli* at 25 µg/ml in liquid or solid media and at 50-200 µg/ml to select Zeocin™-resistant mammalian cells.

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



SgfI (11)
1 GGATCTGCGATCGCTCCGGTGCCCGTCACTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGTTCGCAATTGAACGGTGCCTA
101 GAGAAGTGGCGCGGGTAAACTGGGAAAGTGATGTCGTGTACTGGCTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (246)
PvuII (242)
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCCTTCACGCGCCCGCCCTACCTGAGGCC
301 GCCATCCACGCGGTTGAGTCGCGTTCTGCCGCCTCCCGCTGTGGTGCCTCCTGAATCGCTCCGCGCTAGGTAAGTTTAAAGCTCAGGTCGAGACC
401 GGGCCTTTGTCGGCGCTCCCTTGAGGCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCTGACCCTGCTTGCTCAACTCTACGTCTTTGTTTCGTTT

EcoRI (559) XhoI (570) BglII (586)
AgeI (553) EcoRV (567) NcoI (578)
501 TCTGTTCTGGCCGTTACAGATCCAAGCTGTGACC GGCGCTACCTGAGATCACCGTGAATTCGATATCTCGAGCACCATGGTTAGATCTGTGGAGTGC
1 Val Gl uCys
BbrPI (698)
601 CCACCTTGCCAGCACACCTGTGGCAGGACCTTCAGTCTTCTCTTCCCCCAAACCAAGGACACCCTGATGATCTCCAGAACCCTGAGGTACAGT
4 P roP roCysP roAl aP roP roValAl aGl yP roSer Val P heLeuP heP roP roLysP roLysAspThr LeuMe tI eSer ArgThr P roGl uVal Thr C
701 GCGTGGTGGTGGAGTGGAGCCAGACCCGAGGTCCAGTTCAACTGGTTCAGTGGACGGCATGGAGTGCATAATGCCAAGACAAAGCCCGGGAGGA
37 y sVal Val Val AspVal Ser Hi sGl uAspP roGl uVal Gl nP heAsnTrpTyrVal AspGl yMe tGl uVal Hi sAsnAl aLysThr LysP roArgGl uGl
801 GCAGTTCAACAGCAGCTTCCGTGTGGTCAAGCTCTCACCGTCTGACCAAGGACTGGCTGAACGGCAAGGAGTACAAGTCAAGGTCTCCAACAAAGGC
70 uGl nP heAsnSer Thr P heArgVal Val Ser Val LeuThr Val Val Hi sGl nAspTrpLeuAsnGl yLysGl uTyrLysCysLysVal SerAsnLysGl y
901 CTCCCAGCCCCATCGAGAAAACATCTCCAAAACAAAGGGCAGCCCCGAGAACCACAGGTGTACACCCTGCCCCATCCCGGAGGAGATGACCAAGA
104 P roP roAl aP roI l eGl uLysThr I l eSer LysThr LysGl yGl nP roArgGl uP roGl nVal TyrThr LeuP roP roSer ArgGl uGl uMe tThr LysA
1001 ACCAGGTGAGCTGACCTGCCTGGTCAAAGGCTTACCCAGCGACATCCCGTGGAGTGGGAGAGCAATGGGCAGCCGAGAACCACTACAAGACCAC
137 s nGl nVal Ser LeuThr CysLeuVal LysGl yP heTyrP roSer AspI l eAl aVal Gl uT rpGl uSerAsnGl yGl nP roGl uAsnAsnTyrLysThr Th
1101 ACCTCCATGCTGGACTCCGACGGCTCTTCTCTCTACAGCAAGCTCACCGTGGACAAGAGCAGGTGGCAGCAGGGGAACGTCTTCTCATGCTCCGTG
170 r P roP roMe tLeuAspSer AspGl ySer P heP heLeuTyrSer LysLeuThr Val AspLysSer ArgT rpGl nGl yAsnVal P heSer CysSer Val
XmnI (1183)
NheI (1272)
1201 ATGCATGAGGCTCTGCACAACCACTACACAGAAGAGCCTCTCCCTGTCTCCGGTAAATGAgTgcacgGCTAGCTGGCCAGACATGATAAGATACAT
204 Me tHi sGl uAl aLeuHi sAsnHi sTyrThr Gl nLysSer LeuSer LeuSer P roGl yLys ●●●
1301 TGATGAGTTTGGACAACCAACTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGTGC

HpaI (1412)
1401 AATAACAAGTTAAACAACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTG

AseI (1509)
XmnI (1510)
1501 GTATGGAATTAATTTCAAATACAGCATAGCAAACTTAACTCCAATCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATC
1601 AGGGGCTGTTGCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTCATGGAGTTTAAAGATATAGTGATTTTCCCAAGGTTTGAAGTCTTTCAT

SwaI (1766)
1701 TTCTTTATGTTTTAAATGCAGTACCTCCACATTCCTTTTTAGTAAAATATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTT
1801 ATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTAAATAGAAATGGACAGCAA
1901 GAAAGCGAGCTTCTAGTTATCTCAGTCTGCTCTGCCACAAGTGACAGCAGTTGCCGGCCGGTCCGCGAGGGCGAACTCCCGCCCCACGGCT
125 ●●●AspGl nGl uGl uAl aVal P heHi sVal CysAsnGl yAl aP roAspArgLeuAl aP heGl uArgGl yTrpP roGl
2001 GCTCGCCGATCTCGGTATGGCCGGCCGGAGGCGTCCCGGAAGTTCGTGGACACGACCTCCGACACTCGGGTACAGCTCGTCCAGGCGCGCACCA
99 nGl uGl yI l eGl uThr Me tAl aP roGl ySer Al aAspArgP heAsnThr Ser Val Val Gl uSer TrpGl uAl aTyrLeuGl uAspLeuGl yArgVal T rp
2101 CACCCAGGCCAGGTTGTGTCCGGCACCTGCTGTCGACCGCGCTGATGAACAGGGTTCAGTCTCCCGACCACACCGGCGAAGTCTCTCCACG
66 Val TrpAl aLeuThrAsnAspP roVal Val Gl nAspGl nValAl aSer I l eP heLeuThr Val AspAspArgVal Val Gl yAl aP heAspAspGl uVal P
2201 AAGTCCCGGAGAACCAGCCGGTCCGAGTCCGACCTCCGGCAGCTCGCGCGGGTGAACCCGGAACGGCACTGGTCAACTTGGCCATGA
32 heAspArgSer P heGl yLeuArgAspThr T rpP heGl uValAl aGl yAl aVal AspArgAl aThr LeuVal I P roValAl aSer Thr LeuLysAl aMe t
2301 TGGCTCCTCctgtcaggagaggaagagaagaaggttagtacaattgCTATAGTGAGTTGATTATACTATGCAGATATACTATGCCAATGATTAATTGT
AseI (2393)

PstI (2415)
2401 CAAACTAGGGTGCAGgttcatagtgccacttttctgactgcccactctcctgccaccctttccaggcatagacagtcagtgacttacCAAATC

HindIII (2519)
2501 ACAGGAGGGAGAAGCAGAAGCTTGAGACAGACCCGCGGACCGCCAACTGCGAGGGACGTGGCTAGGGCGGCTTCTTTTATGGTGCGCCGCCCTCG

BspEI (2677)
2601 GAGGAGGGCGCTCGGGAGGCTAGCGCCAATCTGCGGTGGCAGGAGGCGGGCCGAAGCCGTGCCTGACCAATCCGGAGCACATAGGAGTCTCAGC

SpeI (2784)
2701 CCCCCGCCCCAAAGCAAGGGAAGTACGCGCCTGTAGCGCCAGCGTGTGTGAAATGGGGCTTGGGGGTTGGGGCCCTGACTAGTCAAACAAACT
2801 CCCATTGACGTCAATGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCCTATCCACGCCATTGATGTAAGTCCAAAACCGCATCATCATGTAATA

2901 GCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTACCAGTCATTGACGTCAA
3001 TAGGGGGCGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGCAGTTTACCCTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGG

PstI (3200)
SdaI (3200)

3101 CGTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGGTCTGTTGGGCGGTGAGCCAGGCGGGCCATTACCCTAAGTTATGTAACGCTGCA

PacI (3207) **BspLU11I (3213)**

3201 GGTTAATTAAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGA

3301 GCATCACA AAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTGGAAGCTCCCTCGTGCGCTCTCCT

3401 GTTCGACCCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGT

3501 AGGTCGTTGCTCCAAGCTGGCTGTGTGCACGAACCCCGTTAGCCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAG

3601 ACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTAC

3701 GGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCG

3801 CTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCA

PacI (3947) SwaI (3955) **NotI (3963)**

3901 GTGGAACGAAAACACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCCAATAAAATATCTTTATTTTCATTACATCTGT

4001 GTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACAAACTAGCAAATAGGCTGTCCCCAGTGCAAGTGC

4101 AGGTGCCAGAACATTTCTCTATCGAA