

# pSELECT-zeo-mcs

Dual expression cassette plasmid for the expression of one gene of interest

Catalog code: psetz-mcs

For research use only

Version 20L01-MM

## PRODUCT INFORMATION

### Content:

- 20 µg of pSELECT-zeo-mcs plasmid provided as lyophilized DNA
- 1 ml of Zeocin™ (100 mg/ml)

### Storage and Stability:

Product is shipped at room temperature. Lyophilized DNA is stable for 12 months when stored at -20°C. Resuspended DNA is stable for 12 months when stored at -20°C. Avoid repeated freeze-thaw cycles. 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

pSELECT plasmids are specifically designed for strong and constitutive expression of a gene of interest in a wide variety of cell lines. They allow the selection of stable transfectants and offer a variety of selectable markers. pSELECT plasmids contain two expression cassettes: the first drives the expression of the gene of interest and the second drives the expression of a large choice of dominant selectable markers for both *E. coli* and mammalian cells. They are both terminating with a strong polyadenylation signal (polyA) that separates the two expression cassettes thus preventing any transcription interference. The late SV40 polyA terminates the transcription of the gene of interest while the human β-globin polyA terminates the transcription of the selectable marker.

## PLASMID FEATURES

### First expression cassette

- **hEF1-HTLV prom** is a composite promoter comprising the Elongation Factor-1α (EF-1α) core promoter<sup>1</sup> and the R segment and part of the U5 sequence (R-U5') of the Human T-Cell Leukemia Virus (HTLV) Type 1 Long Terminal Repeat<sup>2</sup>. 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 the following restriction sites: 5' - Sal I, SgrA I, BamH I, Eco47 III, Nco I, Nhe I - 3'
- Each restriction site is compatible with many other enzymes, increasing the cloning options.
- **SV40 pAn:** the Simian Virus 40 late polyadenylation signal enables efficient cleavage and polyadenylation reactions resulting in high levels of steady-state mRNA<sup>3</sup>.
  - **Ori:** a minimal *E. coli* origin of replication to limit vector size, but with the same activity as the longer Ori.

## Second expression cassette

- **CMV enh/prom:** The human cytomegalovirus immediate-early gene 1 promoter/enhancer was originally isolated from the Towne strain and was found to be stronger than any other viral promoters.
- **EM7** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*.
- **Zeo:** Resistance to Zeocin™ is conferred by the *Sh ble* gene from *Streptoalloteichus hindustanus*. The *Sh ble* gene is driven by the CMV enhancer/promoter in tandem with the bacterial EM7 promoter allowing 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 transcription<sup>4</sup>.

## 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

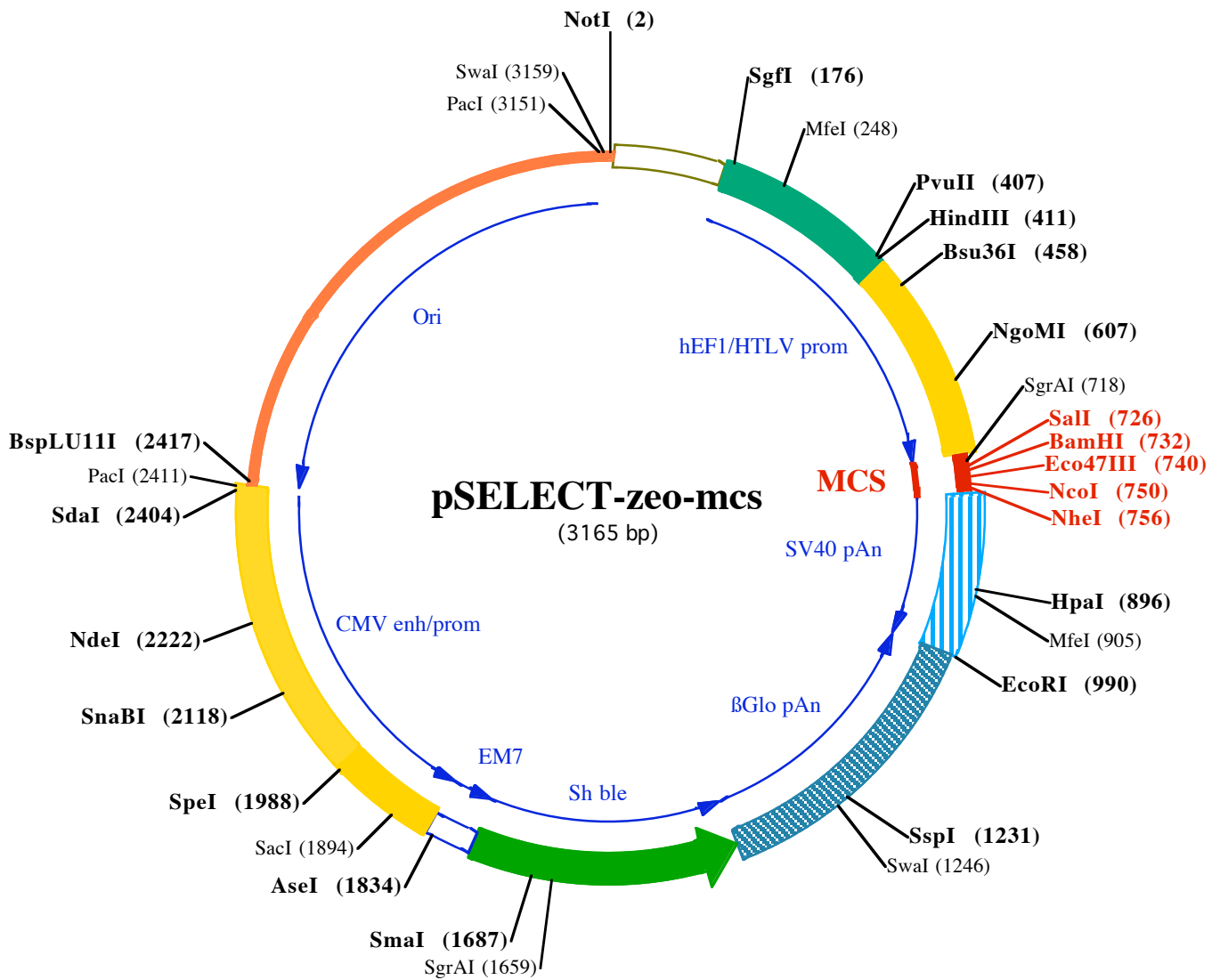
This antibiotic can be used for *E. coli* at 25 µg/ml in liquid or solid media and at 50-200 µg/ml to select Zeocin™-resistant mammalian cells.

### 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](mailto:info@invivogen.com)



90

**NotI (2)**  
1 GCGGCGCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGTTTTTTGTGTGAATCGTAACATAACAGCTCTCCATCAAAACAAAACGAAACA  
**SgfI (176)**  
101 AAACAACTAGCAAATAGGCTGTCCCCAGTGAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGATCTGCGATCGCTCCGGTGCCCGTCAGTGGCCA  
**MfeI (248)**  
201 GAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGTTCGGCAATTGAACGGTGCCTAGAGAAGGTGGCGGGGTAAACTGGGAAAGTGATG  
301 TCGTGTACTGGCTCCGCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCCGTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAG  
**HindIII (411)**  
**PvuII (407)** **Bsu36I (458)**  
401 AACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTTCACGCGCCGCCCTACCTGAGGCGGCCATCCACGCGGTTGAGTCGCGTTCTGCCGCTT  
501 CCGCCTGTGGTGCCTCCTGAAGTGCCTCCGCGTCTAGTTAAGTTTAAAGTCAAGTGCAGACGGGCTTTGTCCGCGCTCCCTTGGAGCTACCTA  
**NgoMI (607)**  
601 GACTCAGCCGGCTCTCCACGCTTTCCTGACCCCTGCTTCTCAACTCTACGTCCTTTGTTTCTGTTTCTGCGCCGTTACAGATCCAAGCTGTGACC  
**BamHI (732)** **NheI (756)**  
**SgrAI (718)** **SalI (726)** **Eco47III (740)** **NcoI (750)**  
701 GCGGCTACCTGAGATCAcggcgtgtcgacggatccagcgtctcgagCCATGGCTAGCTGGCAGACATGATAAGATACATTGATGAGTTTGGACAA  
**HpaI (896)**  
801 ACCACAAC TAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACA  
**MfeI (905)** **EcoRI (990)**  
901 ACAACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGGTATGGAAATCTAAAAA  
1001 TACAGCATAGCAAACTTTAACCTCAAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGC  
1101 ATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCAAAGTTTGAAGTACTCTTCTTTCTTTATGTTTTAAATGCA  
**SspI (1231)** **SwaI (1246)**  
1201 CTGACCTCCACATTCCTTTTTTAGTAAAATATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAAATCCAGATGC  
1301 TCAAGGCCCTTCATAATATCCCCAGTTTGTAGTGTGGACTTAGGGAACAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTA  
1401 TCCTCAGTCTGCTCCTCTGCCACAAAGTGCACGAGTTGCCGGCGGGTTCGCGCAGGGCGAACTCCCGCCACCGGCTGCTCGCGATCTCGGTCATG  
126 GI y ••• AspGI nGI uGI uAl aVal PheHi sVal CysAsnGI yAl aP roAspArgLeuAl aPheGI uArgGI yTrpP roGI nGI uGI yI l eGI uThr Me tA  
1501 GCCGGCCCGGAGGCGTCCCGAAGTTCTGTGGACACGACCTCCGACCACTCGCGGTACAGCTCGTCCAGGCGCGCACCCACCCAGGCCAGGGTGTGT  
92 l aP roGI ySer Al aAspArgPheAsnThr Ser Val Val GI uSer TrpGI uAl aTyrLeuGI uAspLeuGI yArgVal l TrpVal l TrpAl aLeuThrAsnAs  
**SgrAI (1659)** **SmaI (1687)**  
1601 CCGGCACCACTGGTCTGGACCGCGCTGATGAACAGGGTCACTGCTCCCGGACACACCGGCGAAGTCTCTCCACGAAGTCCCGGGAAGACCCGAG  
59 pP roVal Val GI nAspGI nVal Al aSer l l ePheLeuThr Val AspAspArgVal Val GI yAl aPheAspAspGI uVal l PheAspArgSer PheGI yLeu  
1701 CCGGTCGGTCCAGAACTCCGGCAGCTCGCGCGGGTGGACACCGGAACGGCACTGGTCAACTTGGCCATGATGGCCCTCTATAGTGAGTC  
26 l ArgAspThr TrpPheGI uVal Al aGI yAl aVal AspArgAl aThr LeuVal l ProVal Al aSer Thr LeuLysAl aMe t  
**AseI (1834)** **SacI (1894)**  
1801 GTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCACCTAACAGGCTCTGCTT  
**SpeI (1988)**  
1901 ATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAACA  
2001 AACTCCATTGACGCTCAATGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCAGCCATTGATGTACTGCCAAAACCGCATCATCATGGT  
**SnaBI (2118)**  
2101 AATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGTGATGACTGGGCATAATGCCAGGCGGGCCATTACCGTCATTGACC  
**NdeI (2222)**  
2201 TCAATAGGGGGCTACTTGGCATATGATACACTTGTGACTGCCAAGTGGCGAGTTTACCCTAAATACTCCACCCATTGACGTCAATGAAAGTCCCTA  
**SdaI (2404)**  
2301 TTGGCGTTACTATGGGAACATACGTCATTATTGACGCTCAATGGGCGGGGCTGTTGGCGGTGACCCAGGCGGGCCATTACCGTAAGTTATGTAACGCC  
**PaeI (2411)** **BspLU11I (2417)**  
2401 TGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTG  
2501 ACGAGCATCACAAAATCGACGCTCAAGTCAAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCTTTCCCGTGGAAAGTCCCTCGTGGCTC  
2601 TCCTGTCCGACCTCGCCCTTACCGGATACCTGTCCGCTTTCTCCCTTCGGAAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCCG  
2701 GTGTAGGTCGTTGCTCCAAGTGGGCTGTGTGCACGAACCCCGTTCAGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGG  
2801 TAAGACACGACTTATCGCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAA  
2901 CTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAAACAAAC  
3001 ACCGCTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGCTGACG  
**PaeI (3151)** **SwaI (3159)**  
3101 CTCAGTGGAACTCACGTTAAGGATTTTGGTCATGGCTAATTAACATTTAAATCA