

pZERO-mcs

A plasmid containing a multiple cloning site and the puromycin resistance gene

Catalog # pzero-mcs

For research use only

Version # 10C15-MM

PRODUCT INFORMATION

Content:

- 1 disk of lyophilized GT110 bacteria transformed with pZERO-mcs.
- GT110 genotype is: *F⁻, mcrA, Δ(mrr-hsdRMS-mcrBC), Ø80lacΔM15, ΔlacX74, recA1, endA1*.
- 4 pouches of *E. coli* Fast-Media® Puro (2 TB and 2 agar).

Storage and stability:

- Products are shipped at room temperature.
- Transformed bacteria should be stored at -20°C and are stable up to 1 year.
- Store *E. coli* Fast-Media® Puro at room temperature. Fast-Media® pouches are stable 18 months when stored properly.

Quality control:

- Plasmid construct has been confirmed by restriction analysis and sequencing.
- Bacteria have been lyophilized, and their viability upon resuspension has been verified.

GENERAL PRODUCT USE

pZERO-mcs is a ready-made expression vector containing the Puromycin resistance gene, the hybrid EF1α/HTLV promoter and a multiple cloning site. **pZERO-mcs** may be used for:

Cloning in a gene of interest. Six unique restriction sites comprise the MCS facilitating cloning of genes. Cloned genes will be under the control of the EF1α/HTLV promoter.

As an “empty” control vector. Since pZERO-mcs does not contain a gene of interest, it can be used in conjunction with other vectors of the pZERO family to serve as an experimental control.

pZERO-mcs is selectable with the potent antibiotic puromycin in both *E. coli* and mammalian cells.

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 the following restriction sites: 5' - Sal I, BamH I, Eco47 III, Nco I, Nhe I, Msc 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 mRNA³.
- **ori:** a minimal *E. coli* origin of replication to limit vector size, but with the same activity as the longer Ori.
- **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*.

- **Puro:** Resistance to Puromycin is conferred by the *Pac* gene from *Streptomyces* which encodes a N-acetyl-transferase. The *Pac* gene is driven by the CMV enhancer/promoter in tandem with the bacterial EM7 promoter allowing 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 transcription⁴.

METHODS

Growth of pZERO-mcs-transformed bacteria:

Use sterile conditions to do the following:

- 1- Resuspend the lyophilized *E. coli* by adding 1 ml of LB medium in the tube containing the disk. Let sit for 5 minutes. Mix gently by inverting the tube several times.
- 2- Streak bacteria taken from this suspension on a puromycin LB agar plate prepared with the *E. coli* Fast-Media® Puro Agar provided (see below).
- 3- Place the plate in an incubator at 37°C overnight.
- 4- Isolate a single colony and grow the bacteria in TB supplemented with puromycin using the Fast-Media® Puro TB provided (see below).
- 5- Extract the pZERO-mcs plasmid DNA using the method of your choice.

Note: For long-term storage of the pZERO-mcs-transformed bacteria, prepare a 20% glycerol stock of the bacteria grown in the overnight liquid culture and freeze at -80°C.

Selection of bacteria with *E. coli* Fast-Media

Fast-Media® is a **fast and convenient** way to prepare liquid and solid media for bacterial culture by using only a microwave. Fast-Media® is a TB (liquid) or LB (solid) based medium that already contains the antibiotic.

Fast-Media® Puro can be ordered separately (#fas-pr-l (liquid), #fas-pr-s (solid)).

Method:

- 1- Pour the contents of a Fast-Media® pouch into a clean borosilicate glass bottle or flask.
- 2- Add 200 ml of distilled water to the flask
- 3- Heat in a microwave on MEDIUM power setting (about 400Watts), until bubbles start appearing (approximately 3 minutes). **Do not heat a closed container. Do not autoclave Fast-Media®.**
- 4- Swirl gently to mix the preparation. **Be careful, the bottle and media are hot, use heatproof pads or gloves and care when handling.**
- 5- Reheat the media for 30 seconds and gently swirl again. Repeat as necessary to completely dissolve the powder into solution. But be careful to avoid overboiling and volume loss.
- 6- Let agar medium cool to 45°C before pouring plates. Let liquid media cool to 37°C before seeding bacteria.

Note: Do not reheat solidified Fast-Media® as the antibiotic will be permanently destroyed by the procedure.

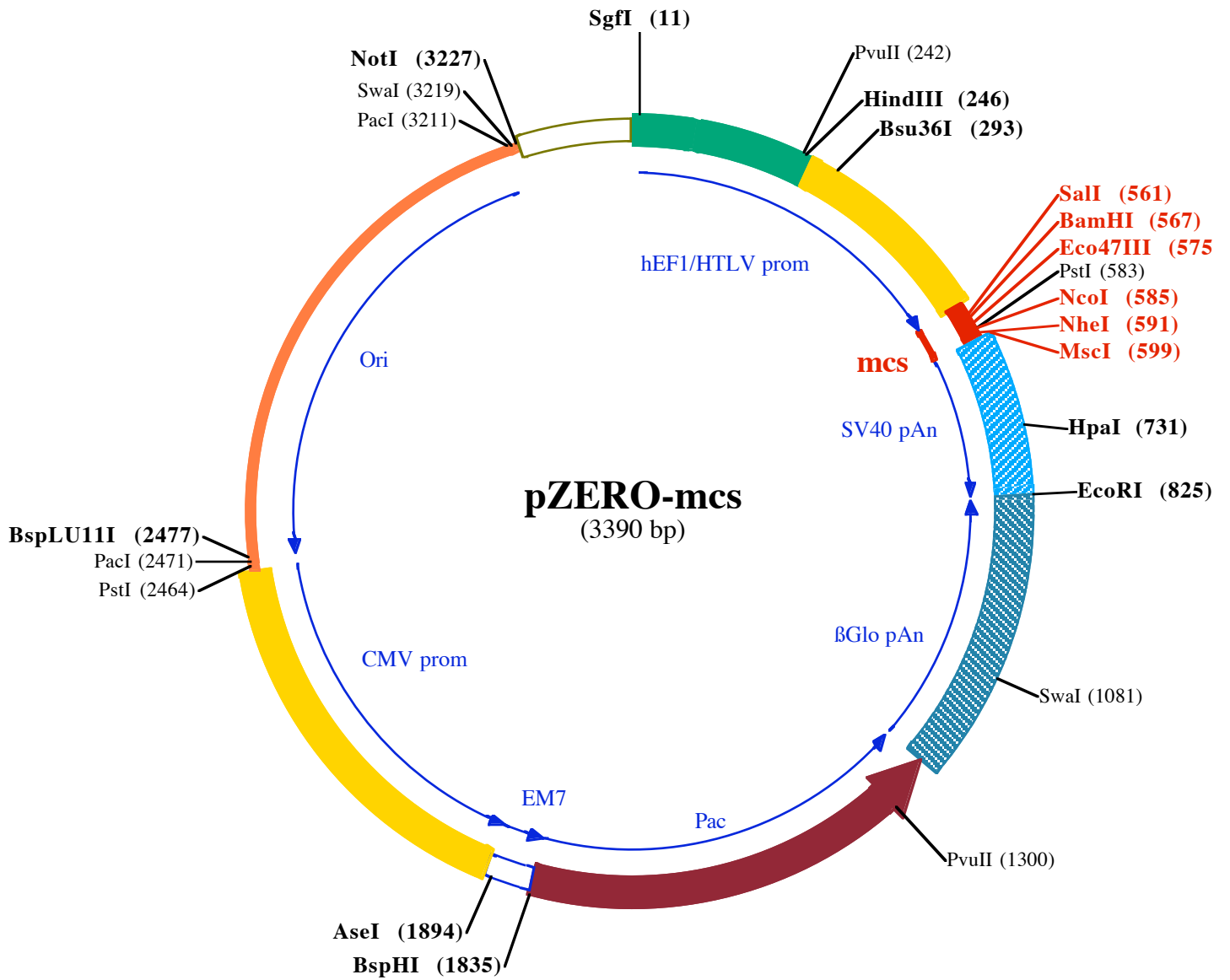
References:

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

Toll free (US): 888-457-5873
Outside US: (+1) 858-457-5873
Europe: +33 562-71-69-39
E-mail: info@invivogen.com
Website: www.invivogen.com


3950 Sorrento Valley Blvd. Suite 100
San Diego, CA 92121 - USA



90

SgfI (11)

1 GGATCTGCGATCGCTCCGGTCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA

101 GAGAAGGTGGCGGGGTAACCTGGAAAGTGATGTCGTGACTGGCTCCGCTTTTCCGAGGGTGGGGAGAACCCTATATAAGTCAGTAGTCGCC

HindIII (246) **PvuII (242)** **Bsu36I (293)**

201 GTGAACGTTCTTTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACGCGCCCGCCCTACCTGAGGCC

301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCTGAACTCGCTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

401 GGGCCTTTGTCGGCGCTCCCTTGGAGCTACCTAGACTCAGCCGGCTCCACGCTTTGCTGACCTGCTTCTCAACTCTACGCTTTTGTTCGTTT

BamHI (567) **PstI (583)** **NheI (591)**
SalI (561) **Eco47III (575)** **NcoI (585)** **MseI (599)**

501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCCCTACCTGAGATCACCGGCGTGTGACGGATCCAGCGCTCTGCAGCCATGGGCTAGCTGGC

601 CAGACATGATAAGATACATTGATGAGTTGGACAAACCAACTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATT

HpaI (731)

701 TGTAACCATTATAAGCTGCAATAACAAGTTAACAACAACAATTGCATTCTTTTATGTTTCAGGTTTCAGGGGAGGTGTGGAGGTTTTTAAAGCAAG

EcoRI (825)

801 TAAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAA

901 TAAGGCATAGGCATCAGGGGCTGTTGCAATGTGCATTAGCTGTTTGACGCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCAAGGTTT

SwaI (1081)

1001 GAACTAGCTCTTCATTTCTTTATGTTTTAAATGCACCTGACCTCCACATTCCCTTTTATGATAAATATTCAGAAATAATTTAAATACATCATTGCAATGA

1101 AAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATA

PvuII (1300)

1201 GAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTCAGGTTTAAAGCTCAGGCTTCTTGTGCATGCACCAAGTCTTGGCCCTTCTGGAACCTCAACATCAG

1301 CTGTCACAGTGAATCCAGTCTTTCATAAAAAGGCAGGTTTCTGGGAGCAGAAGTTCCAGAAAGGCAGGAACCTCAGCCCTTTCAGCAGCTCAACTCC

1794 aThr Val Thr PheGlyLeuArgGlyTyrPheProLeuAsnArgProAlaSer Thr GlyLeuPheAlaProVal GlyAlaArgGlyAlaAlaGlyVal Gly

1401 AGGCAGAACACAGCAGATCCAGACCTTCCCTGGTGGTCAGGGCTCACTCAACAGTTGCCAGAAACCAAGCTGGCTCTTTGGCTGTGTGGTGC

1464 ProLeuVal ValAlaSer GlyLeuGlyLysGlyGlnHisAspProSer Val GlyVal ThrAlaLeuPheTrpAlaProGlyLysProArgHisProAla

1501 AGCAGACCTCCATTTGTTGTTGCTGCCAGCCTGCTCCAGAGAGCTCAGCCATTCTGGTCCAATTCAGCAAAACAGCACCAGCTTCAACAGACT

1124 euLeuGlyGlyuMetGlnGlnAlaAlaLeuArgSer GlySerLeuGlyAlaMetArgProGlyIleGlyAlaPheValAlaGlyAlaGlyValSerGly

1601 CAGTGTTGTCAAACCTGCAACAGCAGCTCCATCATCTGCAACCAACTTTCCAATGTCCAGTCCCACTCTGGTGAGGAAGATTTCTGCAGTTCTGT

794 uProThr Thr TrpValAlaValAlaAlaGlyAspAspAlaVal TrpVal LysGlyIleAspLeuGlyValIleArgThrLeuPheLeuGlyGlnLeuGlyuThr

1701 CACCCTCAATGTGCCTGTGAGGTCAGGTCAGCTGTGCTTGTTCAGGGTAGTCTGCAAAAGCAGCAGCCAGTGTCTCACAGCTCTTGAACATCATCT

464 ValArgGlyuIleHisArgAspProAspValThrHisArgThrAlaProTyrAspAlaPheAlaAlaAlaLeuThrArgValAlaArgProValAspAspAla

BspHI (1835) **AseI (1894)**

1801 CTGGTTGCCAGCCTCACTGTGGTTGTAAGTCACTGATGATGGCCCTCTATAGTGTGCTATTATACTATGCCGATATACTATGCCGATGATTAATTG

124 rGThrAlaLeuArgValThrProLysTyrGlyuThrMet

1901 TCAAACAGCGTGGATGGCTCTCCAGCTTATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCTACCGCCATTGGC

2001 TCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAACAAACTCCCATGACGCTCAATGGGGTGGAGACTTGGAAATCC

2101 CCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGA

2201 AAGTCCCATAAAGTCTACTGTTGGCATAATGCCAGCGGGCCATTTACCCTCATTGACGCTCAATAGGGGGCTACTTGGCATATGATACACTTGTATGTA

2301 CTGCCAAGTGGGAGTTTACCCTAAATACTCCACCCATTGACGCTCAATGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCATTATTGACGTCAA

PacI (2471) **PstI (2464)** **BspLU11I (2477)**

2401 TGGGCGGGGTCGTTGGCGGTCAGCCAGCGGGCCATTTACCCTAAGTTATGTAACCCCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAA

2501 AGGCCAGGAACCGTAAAAAGCCGCTGTGGCGTTTTCCATAGGCTCCGCCCCCTGACGAGCATCAAAAAATCGACGCTCAAGTCAGAGGTGGCG

2601 AAACCCGACAGGACTATAAGATACCAGGCTTTCCCTCGAAGCTCCCTCGTGCCTCTCTGTTCCGACCCTGCCGCTTACCGGATACTGTCCGCC

2701 TTTCTCCTTCGGGAAGCGTGGCGTTTCTCATAGCTCACGCTGATAGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAAC

2801 CCCCCGTTACGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAA

2901 CAGGATTAGCAGAGCGAGGTATGTAGCGGTGTACAGAGTCTTGAAGTGGTGGCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCT

3001 CTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAAACAAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGAAGCAGCAGA

3101 TTACGCGCAGAAAAAAGGATCTCAAGAAGATCTTTGATCTTTTACGGGTCTGACGCTCAGTGAACGAAACTCACGTTAAGGGATTTTGGTCAT

PacI (3211) **SwaI (3219)** **NotI (3227)**

3201 GGCTAGTTAATTAACATTTAAATCAGCGCCGCAATAAAATATCTTTATTTTATTACATCTGTGTGGTTTTTTGTGTAATCGTAACTAACATACG

3301 CTCTCCATCAAAACAAACGAAACAAACAAACTAGCAAAATAGGCTGTCCCGAGTGAAGTGCAGGTGCCAGAACATTTCTCTATCGAA