

# pUNO3-mcs

Expression vector containing a multiple cloning site and the hygromycin resistance gene

Catalog code: puno3-mcs

For research use only

Version 19F21-MM

## PRODUCT INFORMATION

### Contents:

- 20 µg of lyophilized plasmid DNA
- 1 ml Hygromycin B Gold at 100 mg/ml

### Storage and Stability:

- Product is shipped at room temperature.
- Upon receipt, store lyophilized DNA at -20°C.
- Resuspended DNA should be stored at -20°C.
- Store Hygromycin B Gold at 4°C or -20°C. The expiry date is specified on the product label.

### Quality control:

- Plasmid construct has been confirmed by restriction analysis and full-length ORF sequencing.
- Plasmid DNA was purified by ion exchange chromatography.

## GENERAL PRODUCT USE

- **pUNO3-mcs** is a ready-made expression vector containing the hygromycin-resistance gene (*hph*) resistance gene, the hybrid EF1α/HTLV promoter and a multiple cloning site.
- **pUNO3-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.** pUNO3-mcs plasmids were designed to serve as experimental control vectors for the pUNO3 plasmid family.

## 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 water. Store resuspended plasmid at -20°C.

### Plasmid amplification and cloning:

Plasmid amplification and cloning can be performed in *E. coli* GT116 or other commonly used laboratory *E. coli* strains, such as DH5α.

### Hygromycin B usage:

This antibiotic can be used for *E. coli* at 50-100 µg/ml in liquid or solid media and at 50-500 µg/ml to select Hygromycin-resistant mammalian cells.

## PLASMID FEATURES

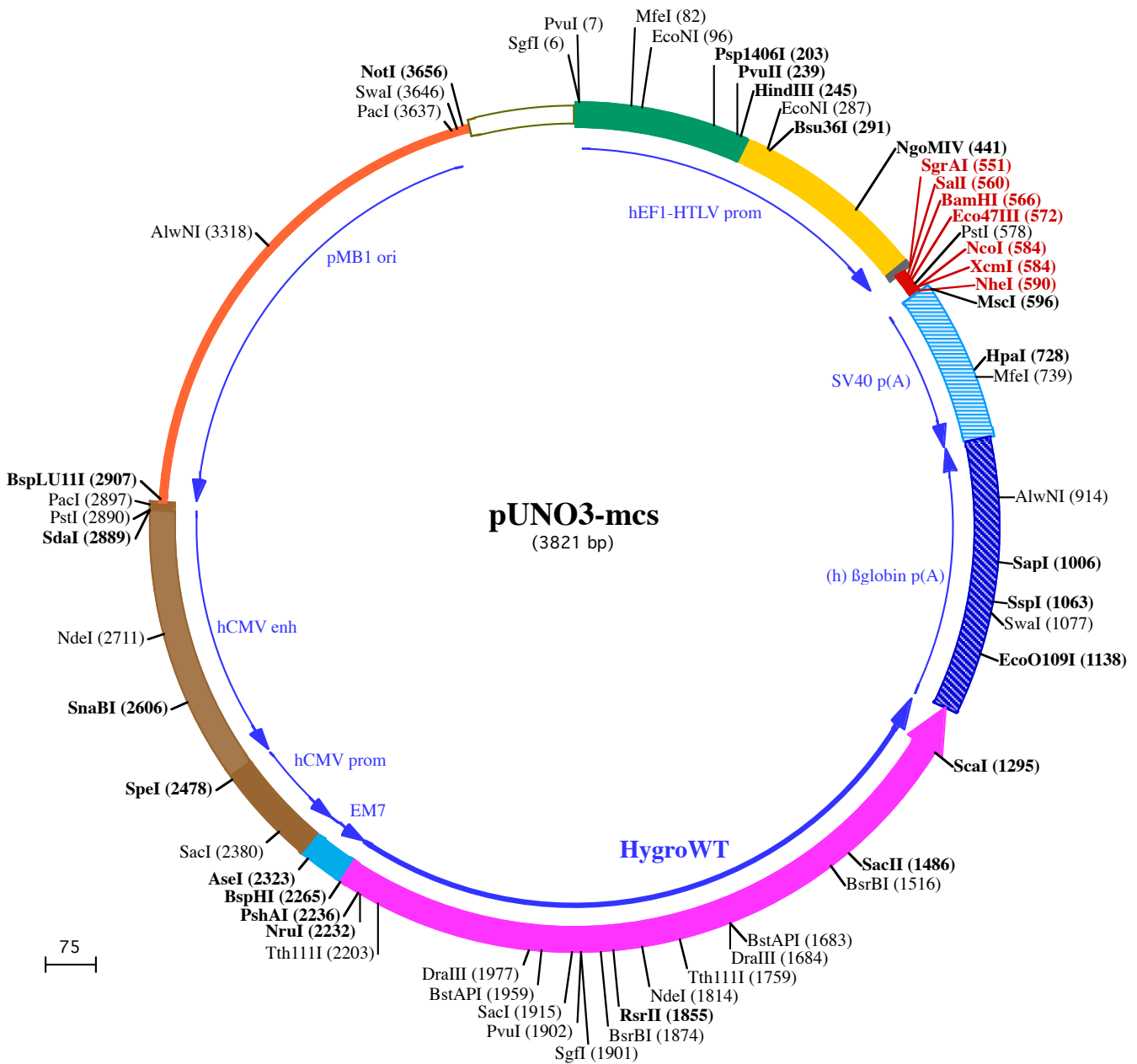
- **EF-1α / HTLV hybrid promoter** is a composite promoter comprised of the Elongation Factor-1α (EF-1α) core promoter<sup>1</sup> and the 5' untranslated region of the Human T-Cell Leukemia Virus (HTLV). EF-1α utilizes a type 2 promoter that encodes for a «house keeping» gene. It is expressed at high levels in all cell cycles and lower levels during G0 phase. The promoter is also non-tissue specific; it is highly expressed in all cell types. The R segment and part of the U5 sequence (R-U5') of the HTLV Type 1 Long Terminal Repeat<sup>2</sup> has been coupled to the EF-1α promoter to enhance stability of DNA and RNA. This modification not only increases steady state transcription, but also significantly increases translation efficiency possibly through mRNA stabilization.
- **MCS:** The multiple cloning site contains the following restriction sites: 5' - SgrAI, Sall, BamHI, Eco47III, NcoI, NheI - 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>.
- **pMB1 ori** is a minimal *E. coli* origin of replication to limit vector size, but with the same activity as the longer Ori.
- **CMV promoter & enhancer** drives the expression of the hygromycin resistance in mammalian cells.
- **Hygro (hygromycin B resistance gene):** Resistance to hygromycin B is conferred by the *hph* gene from *E. coli* which encodes a phosphotransferase. The *hph* gene is driven by the CMV promoter/enhancer in tandem with the bacterial EM7 promoter. Therefore, hygromycin B can be used to select stable mammalian cells transfectants and *E. coli* transformants.
- **Human beta-Globin polyA** is a strong polyadenylation (pAn) signal placed downstream of *hph*. The use of beta-globin pAn minimizes interference<sup>4</sup> and possible recombination events with the SV40 polyadenylation signal.

1. Kim DW. *et al.*, 1990. Use of the human elongation factor 1α promoter as a versatile and efficient expression system. *Gene* 91(2):217-23. 2. Takebe Y. *et al.*, 1988. SR alpha promoter: an efficient and versatile mammalian cDNA expression system composed of the simian virus 40 early promoter and the R-U5 segment of human T-cell leukemia virus type 1 long terminal repeat. *Mol Cell Biol.* 8(1):466-72. 3. Carswell S. & Alwine JC., 1989. Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. *Mol Cell Biol.* 9(10):4248-58. 4. Yu J. & Russell JE., 2001. Structural and functional analysis of an mRNP complex that mediates the high stability of human β-globin mRNA. *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)





PvuI (7) SgfI (6) MfeI (82) EcoNI (96)  
 1 GGATCTGCATCGCTCCGGTGCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA  
 101 GAGAAAGTGGCGCGGGTAAACTGGAAAAGTGTGCTGTACTGGCTCCGCTTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC  
 HindIII (245) Bsu36I (291)  
 Psp1406I (203) PvuII (239) EcoNI (287)  
 201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTCACGCGCCGCGCCCTACCTGAGGCC  
 301 GCCATCCACGCCGGTTGAGTCGCGTTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAACTGCGTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC  
 NgoMIV (441)  
 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTCCTGACCTGCTTCTCAACTCTACGCTTTTGTTCGTTT  
 BamHI (566) PstI (578) XcmI (584) NheI (590)  
 SgrAI (551) Sall (560) Eco47III (572) NcoI (584) MscI (596)  
 501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACTCTGAGATCACCGGCGTGTGACGGATCCAGCGCTCTGCAGCCATGGGCTAGCTGGC  
 601 CAGACATGATAAGATACATTGATGAGTTTGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATT  
 HpaI (728) MfeI (739)  
 701 TGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCATTTTATGTTTCAGGTTCAAGGGGAGGTGTGGGAGTTTTTAAAGCAAG  
 801 TAAACCTCTACAAATGTGGTATGGAATTCTAAATACAGCATAGCAAACCTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAA  
 AlwNI (914)  
 901 TAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGATTAGCTGTTGCAGCCTCACCTTCTTCATGGAGTTAAGATATAGTGATTTTTCCAAGGTTT  
 SapI (1006) SspI (1063) SmaI (1077)  
 1001 GAACTAGCTCTTCATTTCTTTATGTTTTAAATGCAGTACCTCCACATTCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGA  
 EcoO109I (1138)  
 1101 AAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAGGACCTTTAATA  
 ScaI (1295)  
 1201 GAAATTGGACAGCAAGAAAGCGAGCTTCTAGCGAATTCTCGACTATTCTTTGCCCTCGGACGAGTGTGGGGCGTCGGTTTCCACTATCGGCGAGTAC  
 342 E K A R P R T S P R R N G S D A L V  
 1301 TTCTACACAGCCATCGGTCCAGACGGCCGCTTCTCGGGCGATTGTGTACGCCGACAGTCCCGCTCCGGATCGGACGATTGCGTCGCATCGACC  
 323 E V C G D T W V A A S R R A I Q T R G V T G A G S R V I A D C R G  
 SacII (1486)  
 1401 TGCGCCAAGCTGCATCATCGAATTGCCGTCAACCAAGCTCTGATAGAGTTGGTCAAGACCAATGCGGAGCATATACGCCCGGAGCCGCGGATCCTG  
 289 Q A W A A D D F N G D V L S Q Y L Q D L G I R L M Y A R L R P S G A  
 BsrBI (1516)  
 1501 CAAGCTCCGGATGCCTCCGCTCGAAGTAGCGCTGTGCTGCTCATAACAAGCAACCAGGCCTCCAGAAGAAGATGTTGGCGACCTCGATTGGGAATC  
 256 L E P H R R E F Y R T Q Q E M C A L W P R W F F I N A V E Y Q S D  
 DraIII (1684) BstAPI (1683)  
 1601 CCCGAACATCGCCTCGCTCCAGTCAATGACCGCTGTTATGCGGCCATTGTCGCTCAGGACATTGTTGGAGCCGAAATCCGCGTGCACAGGTGCCGGACT  
 223 G F M A E S W D I V A T I R G N D T L V N N S G F D A H V L H R V  
 Tth111I (1759)  
 1701 TCGGGCAGTCTCGGCCAAAGCATCAGCTCATCGAGAGCCTGCGCGACGGACGACTGACGGTGTCTCCATCACAGTTTGCCAGTGATACACATGGG  
 189 E P C D E A W L M L E D L A Q A V S A S V T D D M V T Q W H Y V H P  
 NdeI (1814) RsrII (1855) BsrBI (1874)  
 1801 GATCAGCAATCGCGCATATGAAATCACGCCATGTAGTGTATTGACCGATTCTTCCGGTCCGAATGGGCCGAACCCGCTCGTCTGGCTAAGATCGGCCG  
 156 D A I A C I F D R W T T Y Q G I G Q P G F P G F G S T Q S L D A A  
 PvuI (1902) SgfI (1901) SacI (1915) BstAPI (1959) DraIII (1977)  
 1901 AGCGATCGCATCCATGAGCTCCGCGACGGGTTGAGAACAGCGGGCAGTTCGGTTTTAGGACAGGCTTTCGACCGTGCACCCCTGTGCACGGCGGGAGATG  
 123 A I A D M L E A V P Q L V A P L E T E P L D Q L T V G Q A R R S I  
 2001 CAATAGTTCAGGCTCTCGCTGAATCCCAATGTCAAGCACTTCGGAATCGGAGCGCGCCGATGCAAAGTCCGATAAACATAACGATCTTTGTAGA  
 89 C Y T L S E S F E G I D L V E P I P L A A S A F H R Y V Y R D K Y F  
 2101 AACCATCGGCGCAGCTATTTACCCGAGGACATATCCACGCCCTCTACATCGAAGCTGAAAGCACGAGATTTCTCGCCCTCCGAGAGCTGCATCAGGTC  
 56 G D A C S N V R L V Y G R G G V D F S F A R S E E G E S L Q M L D  
 PshAI (2236) NruI (2232) BspHI (2265)  
 Tth111I (2203)  
 2201 GGAGACGCTGTGAACTTTTCGATCAGAAATTCGCGACAGACGTGCGGTGAGTTGAGGCTTTTTTTCATGATGGCCCTCTATAGTGAGTCGTATTATAC  
 23 S V S D F K E I L F K A V S T A T L E P K K M

2301 TATGCCGATATACTATGCCGATGATTAATTGTCAAACACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACC  
AseI (2323) SacI (2380)

2401 TCCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCCAT  
SpeI (2478)

2501 TGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGAT

2601 GACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGG  
SnaBI (2606)

2701 GGCGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTA  
NdeI (2711)

2801 CTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTGAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACCGCTGCAGGTTA  
PacI (2897)  
PstI (2890)  
SdaI (2889)

2901 ATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATC  
BspLU11I (2907)

3001 ACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCGTGAAGCTCCCTCGTGCCTCTCCTGTTCC

3101 GACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTC

3201 GTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACAG

3301 ACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTA  
AlwNI (3318)

3401 CACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGT

3501 AGCGGTGGTTTTTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCTTTGATCTTTTACGGGTCTGACGCTCAGTGGA

3601 ACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAATCAGCGGCCGCAATAAAAATATCTTTATTTTCATTACATCTGTGTGT  
PacI (3637) SmaI (3646) NotI (3656)

3701 GGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCCGAGTGCAAGTGCAGGTG

3801 CCAGAACATTTCTATCGAA