

pUNO3-mcs

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

Catalog # puno3-mcs

For research use only

Version # 15J27v40-JC

PRODUCT INFORMATION

Contents:

- 20 µg of pUNO3-mcs provided as lyophilized plasmid DNA
- 4 pouches of *E. coli* Fast-Media® Hygro (2 TB and 2 Agar)

Storage and Stability:

- Product is shipped at room temperature.
- Store lyophilized DNA at -20 °C.
- Resuspended DNA is stable for up to 1 year when stored at -20 °C .
- Store *E. coli* Fast-Media® at room temperature in a dry and cool place. Fast-Media® pouches are stable for 2 years when stored properly.

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.

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

E. coli Fast-Media® Hygro is a fast and convenient way to prepare liquid and solid media for bacterial culture by using only a microwave.

1- Pour the contents of a 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 400 Watts), 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.

PLASMID FEATURES

• **EF-1α/HTLV hybrid promoter** is a composite promoter comprised of the Elongation Factor-1α (EF-1α) core promoter¹ 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 Repeat² 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 mRNA³.

• **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.

• **EM7** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*.

• **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 interference⁴ 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.

RELATED PRODUCTS

Product	Catalog Code
Hygromycin B Gold (solution)	ant-hg-1
ChemiComp GT116	gt116-11
Fast-Media™ Hygro Agar	fas-hg-s
Fast-Media™ Hygro TB	fas-hg-l

TECHNICAL SUPPORT

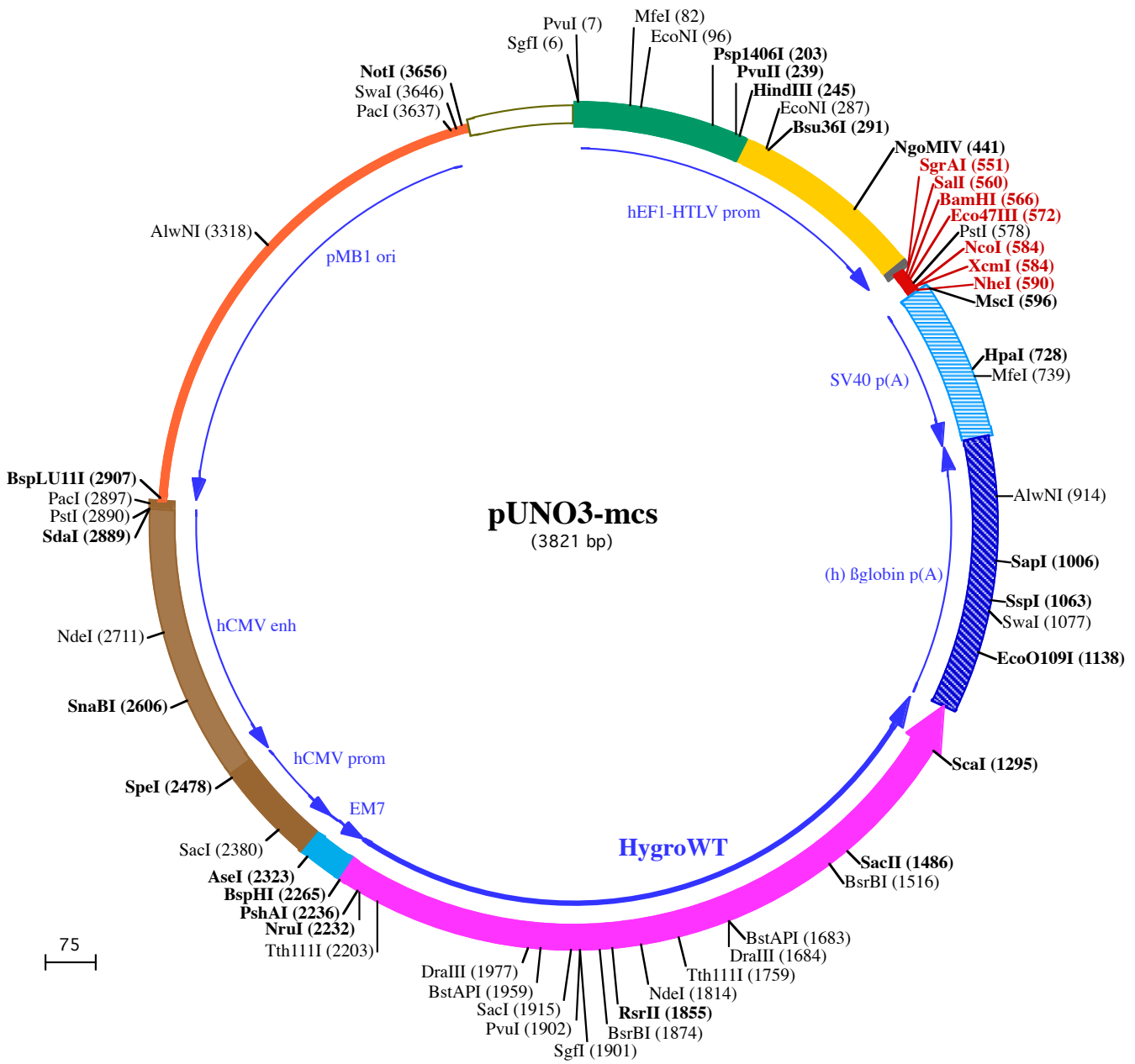
InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3-622-34-80

E-mail: info@invivogen.com



PvuI (7) SgfI (6) MfeI (82) EcoNI (96)
 1 GGATCTGCGATCGCTCCGGTGCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCCGCAATTGAACGGGTGCCTA
 101 GAGAAAGTGGCGCGGGTAAACTGGAAAAGTGTGCTGTACTGGCTCCGCTTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC
 HindIII (245) Bsu36I (291)
 Psp1406I (203) PvuII (239) EcoNI (287)
 201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTCACGCGCCGCCCTACCTGAGGCC
 301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTAAAGCTCAGGTCGAGACC
 NgoMIV (441)
 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCTGACCTGCTTCTCAACTCTACGCTTTTGTTCGTTT
 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 TGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCATTTTATGTTTCAGGTTCCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAG
 801 TAAACCTCTACAAATGTGGTATGGAATTCTAAATACAGCATAGCAAACCTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAA
 AlwNI (914)
 901 TAAGGCATAGGCATCAGGGGCTGTTGCCAATGTCATTAGCTGTTGCAGCCTCACCTTCTTCATGGAGTTAAGATATAGTGATTTTTCCCAAGGTTT
 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 TTCTACACAGCCATCGGTCCAGACGGCCGCGCTTCTCGGGCGATTGTGTACGCCGACAGTCCCGCTCCGGATCGGACGATTGCGTCGCATCGACC
 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 TGCGCCAAGCTGCATCATCGAATTGCCGTCAACCAAGCTCTGATAGAGTTGGTCAAGACCAATGCGGAGCATATACGCCGGAGCCGCGGATCCTG
 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 CAAGTCCGGATGCTCCGCTCGAAGTAGCGCTGTGCTGCTCATAACAAGCAACCAGGCCTCCAGAAGAAGATGTTGGCGACCTCGATTGGGAATC
 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 TCGGGCAGTCTCGGCCAAAGCATCAGCTCATCGAGAGCCTGCGGACGGACGACTGACGGTGTCTCCATCACAGTTTGCCAGTGATACACATGGG
 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 GATCAGCAATCGCGCATATGAAATCACGCCATGTAGTGTATTGACCGATTCTTGGCGTCCGAATGGGCCGAACCCGCTCGTCTGGCTAAGATCGGCCG
 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 AGCGATCGCATCCATGAGCTCCGCGACGGGTTGAGAACAGCGGGCAGTTCGGTTTTAGGCGAGGTTGCAACGTGACACCCTGTGCACGGCGGGGAGATG
 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 CAATAGTTCAGGCTCTCGCTGAATCCCAATGTCAAGCACTTCGGAATCGGGAGCGCGCCGATGCAAAGTCCGATAAACATAACGATCTTTGTAGA
 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 GGAGACGCTGTCGAACTTTTCGATCAGAAATTCGCGACAGACGTCCGGTGAGTTGAGGCTTTTTTTCATGATGGCCCTCTATAGTGAGTCGTATTATAC
 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 TGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTAAGTCCAAAACCGCATCATCATGGTAATAGCGAT

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 CACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCCGCTGGT

3501 AGCGGTGGTTTTTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCTTTGATCTTTTACGGGTCTGACGCTCAGTGGA

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

3701 GGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCCGAGTCAAGTGCAGGTG

3801 CCAGAACATTTCTCTATCGAA