

pUNO2-mcs

A plasmid containing a multiple cloning site and the Zeocin™ resistance gene

Catalog code: puno2-mcs

<https://www.invivogen.com/puno-mcs>

For research use only

Version 21EL19-MM

PRODUCT INFORMATION

Contents

- 20 µg of lyophilized plasmid DNA
- 1 ml of Zeocin™ at 100 mg/ml

Storage and Stability

- Product is shipped at room temperature.
 - Lyophilized DNA should be stored at -20°C.
 - Resuspended DNA should be stored at -20°C and is stable at least for 1 year.
 - Store Zeocin™ 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 open reading frame (ORF) sequencing.
- Plasmid DNA was purified by ion exchange chromatography.

GENERAL PRODUCT USE

pUNO2-mcs is a ready-made expression vector containing the Zeocin™ resistance gene, the hybrid EF1α/HTLV promoter and a multiple cloning site.

pUNO2-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. pUNO2-mcs plasmids were designed to serve as experimental control vectors for the pUNO2 plasmid family that feature a wide choice of native and fusion genes.

METHODS

Plasmid resuspension

Briefly 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.

Zeocin™ usage

Zeocin™ should be used at 25 µg/ml in bacteria and 50-400 µg/ml in mammalian cells. Zeocin™ is supplied as 100 mg/ml solution in HEPES buffer.

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 unique restriction sites:

5' - Sall, BamHI, Eco47III, NcoI, XcmI, 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 Zeocin™ resistance in mammalian cells.

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

• **Sh ble (Zeocin™ resistance gene):** Resistance to Zeocin™ is conferred by the *Sh ble* gene from *Streptoalloteichus hindustanus*. The *Sh ble* gene is driven by the CMV promoter/enhancer in tandem with the bacterial EM7 promoter. Therefore, Zeocin™ 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 *Sh ble*. 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	Description	Cat. Code
Zeocin™	Selection antibiotic	ant-zn-1
ChemiComp GT116	Competent <i>E. coli</i>	gt116-11

TECHNICAL SUPPORT

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

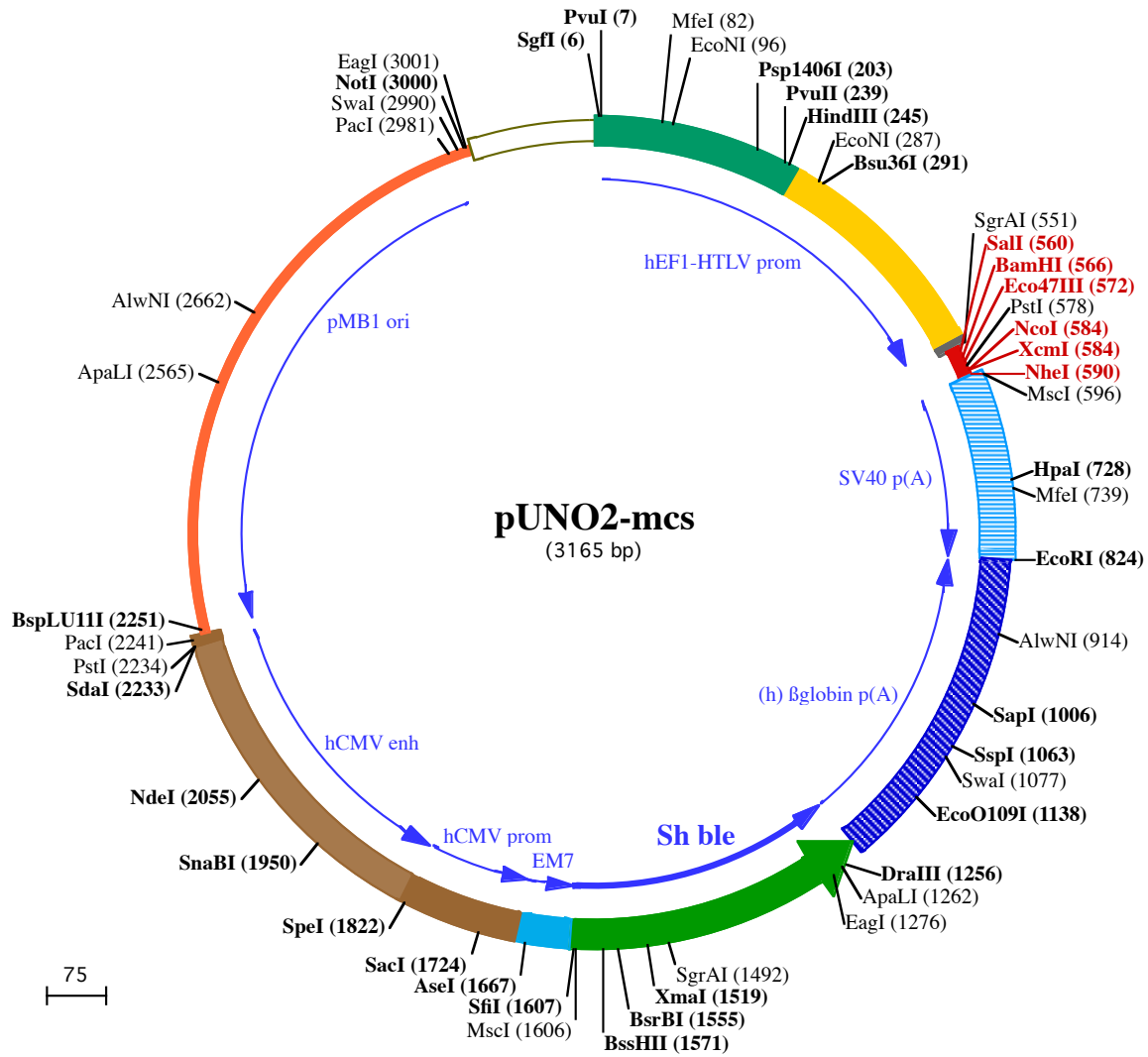
InvivoGen USA (International): +1 (858) 457-5873

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

InvivoGen Asia: +852 3622-3480

E-mail: info@invivogen.com

 **InvivoGen**
www.invivogen.com



PvuI (7)
SgfI (6)
MfeI (82) **EcoNI (96)**

1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA

101 GAGAAGGTGGCGCGGGTAAACTGGAAAAGTATGTCGTGTACTGGCTCCGCCTTTTTCCCGAGGGTGGGGGAGAACCGTATATAAGTGCAGTAGTCGCC

Psp1406I (203)
HindIII (245) **Bsu36I (291)**

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCCTTCACGCGCCCGCCGCTACCTGAGGGCC

301 GCCATCCACGCCGGTTGAGTCGCGTTTCTGCCGCTCCCGCCTGTGGTGCCTCCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTCCTGACCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

XcmI (584)

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATC**BamHI (566)** **PstI (578)** **NheI (590)**
SgrAI (551) **SalI (560)** **Eco4VII (572)** **NcoI (584)** **MscI (596)**

601 CAGACATGATAAGATACATTGATGAGTTTGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATT

HpaI (728) **MfeI (739)**

701 TGTAACCATTATAAGCTGCAATAAAACAAGTTAAACAACAACCTGCATTCTTTATGTTTCAGGTTCCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAG

EcoRI (824)

801 TAAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCAAACCTTAACTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAA

901 TAAGGCATAGGCATCAGGGGCTGTTGCCAATGTCATTAGCTGTTGACGCCTCACCTTCTTCATGGAGTTAAGATATAGTGATTTTTCCCAAGTTTT

AlwNI (914)

1001 GAACTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGA

SapI (1006)
SspI (1063) **Swal (1077)**

1101 AAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATA

EcoO109I (1138)

1201 GAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTCTGCCACAAAGTGACGCAGTTGCCGCGGGGTCCGCGAGGGCGAACT

1301 CCCGCCCCACGGCTGCTCGCCGATCTCGGTATGGCCGGCCCGGAGGCGTCCCGAAAGTTCGTGGACACGACCTCCGACCACTCGCGGTACAGCTCGTC
104 R G W P Q E G I E T M A P G S A D R F N T S V V E S W E A Y L E D

1401 CAGGCCGCGCACCCACCCAGGCCAGGGTGTGTCCGGCACCACTGGTCTGGACCGCGTGATGAACAGGGTCACGTCGTCGCGGACCAACCCGGCGG
71 L G R V W V W A L T N D P V V Q D Q V A S I F L T V D D R V V G A

1501 AAGTCGTCTCCACGAAGTCCCGGGAGAACCCGAGCCGGTCCGATCCAGAACTCGACCGCTCCGGCGACGTCGCGCGCGGTGAGCACCGGAACGGCACTGG
37 F D D E V F D R S F G L R D T W F E V A G A V D R A T L V P V A S T

SfiI (1607)
MscI (1606) **AseI (1667)**

1601 TCAACTTGGCCATGATGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCC
4 L K A M

SacI (1724)

1701 AGCTTATCTGACGGTTCATAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACA

SpeI (1822)

1801 TTTTGGAAAGTCCGTTGATTTACTAGTCAAAAACAACTCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGCTATCCACGC

SnaBI (1950)

1901 CCATTGATGACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCCATAGGTCATGTACTGGG

NdeI (2055)

2001 CATAATGCCAGGCGGGCATTACCCTGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGATGACTGCCAAGTGGGCAGTTTACCCTAA

2101 ATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGAAACATACGTATTGACGTCAATGGGGGGGGTCTTGGCGGTACG

PacI (2241)

PstI (2234)

SdaI (2233) BspLU11I (2251)

2201 CCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGCCGC

2301 GTTGTGGCGTTTTTCCATAGGCTCCGCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATAC

2401 CAGGCGTTTTCCCCTGGAAGCTCCCTCGTGCGCTCTCCTGTTCCGACCTGCCGCTTACCGGATACCTGTCCGCCTTTCTCCCTTCGGGAAGCGTGGCGC

ApaLI (2565)

2501 TTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTTCAGCCGACCGCTGCGC

AlwNI (2662)

2601 CTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTA

2701 GGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAA

2801 AAAGAGTTGGTAGCTCTTGATCCGGCAAACAACCACCGCTGGTAGCGGTGGTTTTTTTGTTCGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCA

PacI (2981) SmaI (2990)

2901 AGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA

EagI (3001)

NotI (3000)

3001 GCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGGTTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACA

3101 AAACAACTAGCAAATAGGCTGTCCCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA