

pMXs-miR-GFP/Puro Retroviral Expression Vector

CATALOG NUMBER: RTV-017

STORAGE: -20°C

QUANTITY AND CONCENTRATION: 10 µg at 0.25 µg/µL in TE

Background

MicroRNAs (miRNAs) are 18–24 nucleotide RNA molecules that regulate the stability or translational efficiency of target mRNAs. These regulatory RNAs function by acting as sequence-specific guides which recruit a large protein complex known as the RNA-induced silencing complex (RISC) to target mRNAs which are subsequently silenced. Diverse functions have been attributed to miRNAs including the regulation of cellular differentiation, proliferation, and apoptosis. Moreover, significant evidence has accumulated implicating a fundamental role for miRNAs in the development of cancer.

miRNAs are initially transcribed as long precursor transcripts known as primary microRNAs (pri-miRNAs). Within these transcripts, the mature miRNA sequences are found in ~60–80 nucleotide hairpin structures. Mature miRNAs are generated from pri-miRNAs by sequential processing (Figure 1). Pri-miRNAs are initially recognized in the nucleus by the microprocessor complex which includes as core components the RNase-III enzyme Drosha and its obligate partner DGCR8. This complex excises the hairpin structure containing the mature miRNA sequence. The liberated hairpins, referred to as precursor miRNAs (pre-miRNAs), are recognized by the nuclear export factor exportin 5 which transports them to the cytoplasm. There, the RNase-III enzyme Dicer performs a second cleavage to generate a double-stranded 18–24 nucleotide RNA molecule. The RISC then associates with this RNA duplex and unwinds it. Generally, only one strand is stably incorporated into the RISC; the other is discarded and rapidly degraded. miRNAs guide the RISC to target messages that are subsequently cleaved or translationally silenced.

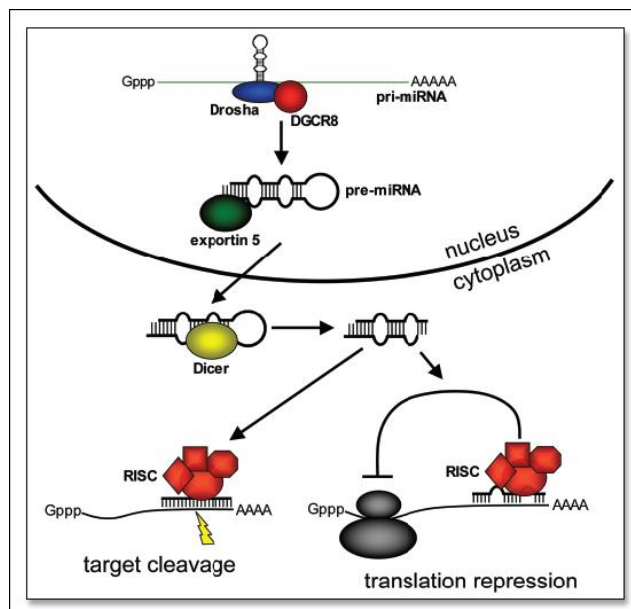


Figure 1. miRNA Biogenesis and function

Retroviruses are efficient tools for delivering heritable genes into the genome of dividing cells. Cell Biolabs' pMXs retroviral vector is based on Moloney murine leukemia virus (MMLV). The vector provides the viral package signal, transcription and processing elements, and MCS for cloning of a target gene. The viral *env* gene, produced by the package cell line, encodes the envelope protein, which determines the viral infectivity range. Transfection into a package cell line produces high-titer, replication-incompetent viruses. In addition to transfer and expression of exogenous genes in mammalian cells, recently, retroviruses have been used to express silencing RNAs (siRNA) to decrease the expression of target genes both *in vitro* and *in vivo*.

Cell Biolabs' pMXs-miR-GFP/Puro Retroviral Expression Vector is designed to clone and express an individual miRNA precursor in its native context while preserving putative hairpin structures to ensure biologically relevant interactions with endogenous processing machinery and regulatory partners, leading to properly cleaved microRNAs. Individual miRNA precursor from any species can be cloned between Xho I sites (Figure 2).

The pMXs-miR-GFP/Puro retroviral expression vector contains the following features:

- **miRNA Processing** – miRNA stem loop precursor in its native context is cloned between Xho I sites. To preserve the putative hairpin structure and proper endogenous processing, miRNA stem loop sequence is flanked by its native intron sequence.
- **EF-1 α Promoter** - ensures a high level of expression in mammalian cells
- **GFP-Puro Fusion Marker** - to monitor cells positive for expression and stable selection with either GFP or puromycin resistance.
- **pUC Origin** - for high copy replication and maintenance of the plasmid in *E. coli*
- **Ampicillin Resistance Gene** - for selection in *E. coli*

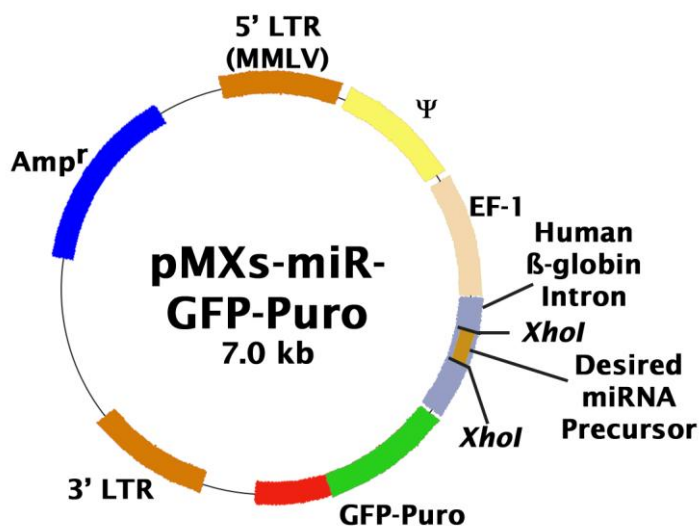
Related Products

1. RV-101: Platinum-E Retroviral Packaging Cell Line, Ecotropic
2. RV-102: Platinum-A Retroviral Packaging Cell Line, Amphotropic
3. RV-103: Platinum-GP Retroviral Packaging Cell Line, Pantropic
4. VPK-120: QuickTiter™ Retrovirus Quantitation Kit
5. RV-200: ViraDuctin™ Retrovirus Transduction Kit

Safety Consideration

Remember that you will be working with samples containing infectious virus. Follow the recommended NIH guidelines for all materials containing BSL-2 organisms. Always wear gloves, use filtered tips and work under a biosafety hood.

pMXs-miR-GFP/Puro Retroviral Expression Vector



MCS: GATTAGTTCTCGAGGATCCGACTGAAGTCGCTAGCTCGAGCTTTTGA
XhoI XhoI

Figure 2: pMXs-miR-GFP/Puro Retroviral Expression Vector (7001 bp, Ampicillin-resistant).

miRNA Precursor Cloning

All of our premade human and mouse miRNA precursor clones in pEP-miR and pEGP-miR vectors are based on the following design, and the resulting overexpression of the mature miRNA is confirmed by Northern blot or Real Time PCR. Here we use human let-7a-2 miRNA as an example:

1. Download desired miRNA stem loop sequence from Sanger's miRNA database:
<http://microrna.sanger.ac.uk/sequences/>

Homo sapiens let-7a-2 stem-loop structure

```
uu  g  u          uagaa ua a
agg gag uag agguuguauaguu    u  c u
||| ||| ||| ||||| ||||| ||| | | c
ucc uuc auc uccgacaugucaa    a  g a
-u  g  c          --uag gg a
```

Homo sapiens let-7a-2 stem-loop sequence

AGGUUGAGGUAGUAGGUUGUAUAGUUUAGAAUUACAUCAAGGGAGAUAACUGUACAGCCUCCUAGCUUUCU

2. Blast search miRNA stem loop sequence: <http://blast.ncbi.nlm.nih.gov/Blast.cgi>

> [ref|NT_033899.7|Hs11_34054](#) D Homo sapiens chromosome 11 genomic contig, reference assembly
Length=38509590

```
Query 1          AGGTTGAGGTAGTAGGTTGTATAGTTTAGAATTACATCAAGGGAGATAACTGTACAGCCT 60
                |||
Sbjct 25579717  AGGTTGAGGTAGTAGGTTGTATAGTTTAGAATTACATCAAGGGAGATAACTGTACAGCCT
25579658

Query 61          CCTAGCTTTCCT 72
                |||
Sbjct 25579657  CCTAGCTTTCCT 25579646
```

3. PCR and Cloning:

- 1) Add 100 base native flank sequence to both upstream and downstream of the miRNA stem loop.

Human let-7a-2 miRNA precursor sequence including the 100 base flank sequences on both ends of the stem loop: let-7a-2 stem-loop sequence is underlined.

```
GCCCAAATAGGTGACAGCACGATGAATCATTATAAGACTAACTTGTAATTTCCCTGCTTAAGAAATG
GTAGTTTTCCAGCCATTGTGACTGCATGCTCCCAGGTTGAGGTAGTAGGTTGTATAGTTTAGAATTA
CATCAAGGGAGATAACTGTACAGCCTCCTAGCTTTCCTTGGGTCTTGCACTAAACAACATGGTGAGA
ACGATCATGATTCCCTCCAGGCCTTTTCTCCCTATGAAAGGTAAGATTGGGTACGATTATTTTATGGT
ATTT
```

- 2) Design PCR primers including Xho I site with four extra bases.

Forward PCR Primer: tcga-ctcgag (XhoI) -21 nt
Reverse PCR Primer: tcga-ctcgag (XhoI) -21 nt

For human let-7a-2 miRNA precursor:

Forward PCR Primer: tcga-ctcgag-gcccaaataaggtagacagcacg
Reverse PCR Primer: tcga-ctcgag-aaataccataaaataatcgta

- 3) PCR the miRNA precursor from genomic DNA and clone into the XhoI sites of the expression vector.

PCR Product of let-7a-2 precursor: let-7a-2 stem-loop sequence is underlined.

```
1 tcgactcgag gcccaaatag gtgacagcac gatgaatcat tataagacta acttgtaatt
61 tccctgctta agaaatggta gttttccagc cattgtgact gcatgctccc aggttgaggt
121 agtaggttgt atagtttaga attacatcaa gggagataac tgtacagcct cctagctttc
181 cttgggtcct gcactaaaca acatgggtgag aacgatcatg attcctccag gcctttttctc
241 cctatgaaaag gtaagattgg gtacgattat tttatgggat ttctcgagtc ga
```

- 4) Validate the insert by DNA sequencing.

Forward Sequencing Primer: TTTGCACCATTCTAAAGAAT
Reverse Sequencing Primer: AAACCTCTTACATCAGTTAC

References

1. microRNA sequences listed in Sanger's miRBase (<http://microrna.sanger.ac.uk/sequences/>).
2. John, B., C. Sander and D. S. Marks (2006) *Methods Mol Biol* **342**: 101-13.
3. Kitamura T., et al., (2003) *Exp. Hematol.* **31**, 1007-1014.

Recent Product Citations

1. Kariba, Y. et al. (2020). Brown adipocyte-derived exosomal miR-132-3p suppress hepatic Srebf1 expression and thereby attenuate expression of lipogenic genes. *Biochem Biophys Res Commun.* S0006-291X(20)31011-1. doi: 10.1016/j.bbrc.2020.05.090.
2. Pollutri, D. et al. (2018). The epigenetically regulated miR-494 associates with stem-cell phenotype and induces sorafenib resistance in hepatocellular carcinoma. *Cell Death Dis.* **9**(1):4. doi: 10.1038/s41419-017-0076-6.
3. Ribocco-Lutkiewicz, M. et al. (2016). MicroRNA expression in amniotic fluid cells. *Fetal Stem Cells in Regenerative Medicine.* doi:10.1007/978-1-4939-3483-6_11.
4. Mansour, M. et al. (2013). The TAL1 complex targets the FBXW7 tumor suppressor by activating miR-223 in human T cell acute lymphoblastic leukemia. *J. Exp. Med.* **210**:1545-1557.

Appendix

Vector Features:

320-614:	MuLv 5' LTR
615-1519:	extended packaging signal (Ψ +)
1520-1851:	Mo-MuLV pol region containing splice acceptor
2049-2444 :	EF-1 promoter
2477-2967:	Human beta-globin inton
2680-2727:	<u>MCS</u>
2995-4317	GFP-Puro Fusion
4384-4975:	MuLv 3' LTR
6028-6888:	Ampicillin Resistance

```
CCCGAAAAGTGCCACCTGCATAATGAAAAGACCCACCTGTAGGTTTGGCAAGCTAGCTTAAGTAACGCCATTTTGCAAGGCATGG
AAAAATACATAACTGAGAATAGAAAAGTTCAGATCAAGGTCAGGAACAGATGGAACAGCTGAATATGGGCCAAACAGGATATCTG
TGGTAAGCAGTTCCTGCCCCGGCTCAGGGCCAAGAACAGATGGAACAGCTGAATATGGGCCAAACAGGATATCTGTGGTAAGCAG
TTCCTGCCCCGGCTCAGGGCCAAGAACAGATGGTCCCCAGATGCGGTCCAGCCCTCAGCAGTTTCTAGAGAACCATCAGATGTTTT
CCAGGGTGCCCCAAGGACCTGAAATGACCCTGTGCCTTATTTGAACTAACCAATCAGTTCGCTTCTCGCTTCTGTTCGCGCGCTT
CTGCTCCCCGAGCTCAATAAAAAGAGCCCAACCCCTCACTCGGCGCGCCAGTTCCTCCGATTGACTGAGTCGCCGGGTACCCGT
GTATCCAATAAACCTCTTGCAGTTGCATCCGACTTGTGGTCTCGCTGTTCCCTGGGAGGGTCTCCTCTGAGTGATTGACTACCC
GTCAGCGGGGGTCTTTCATTTGGGGGCTCGTCCGGGATCGGGAGACCCCTGCCAGGGACCACCGACCCACCACCGGGAGGTAAG
CTGGCCAGCAACTTATCTGTGTCTGTCCGATTGTCTAGTGTCTATGACTGATTTTATGCGCCTGCGTCCGTTACTAGTTAGCTAAC
TAGCTCTGTATCTGGCGGACCCGTGGTGGAACTGACGAGTTTCGGAACACCCGGCCGCAACCTGGGAGACGTCCAGGGACTTCG
GGGCGCTTTTGTGGCCGACCTGAGTCCAAAAATCCCGATCGTTTTGGACTCTTTGGTGCACCCCTAATAGGAGGGATATG
TGGTCTGGTAGGAGACGAGAACCTAAAACAGTTCCCGCCTCCGCTGAATTTTGGCTTTCGGTTTGGGACCGAAGCCGCGCCG
GCGTCTTGTCTGCTGCAGCATCGTTCTGTGTTGTCTCTGTCTGACTGTGTTTTCTGTATTTGTCTGAAAATTAGGGCCAGACTGTT
ACCACTCCCTTAAGTTTACCTTAGGTCACTGGAAAAGATGTCGAGCGGATCGCTCACAACCAGTCGGTAGATGTCAAGAAGAGAC
GTTGGGTTACCTTCTGCTCTGCAGAAATGGCCAACCTTTAACGTCGGATGGCCGCGAGACGGCACCTTTAACCGAGACCTCATCAC
CCAGGTTAAGATCAAGGCTTTTTCACCTGGCCCGCATGGACACCCAGACCAGGTCCCTACATCGTGACCTGGGAAGCCTTGGCT
```

TTTGACCCCTCCCTGGGTCAAGCCCTTTGTACACCCTAAGCCTCCGCTCCTCTTCTCCATCCGCCCCGTCTCTCCCTTG
AACCTCCTCGTTGACCCCGCTCGATCCTCCCTTTATCCAGCCCTCACTCCTTCTTAGGCGCCCCATATGGCCATATGAGAT
CTTATATGGGGCACCCCGCCCTTGTAACTTCCCTGACCTGACATGACAAGAGTTACTAACAGCCCTCTCTCCAAGCTCAC
TTACAGGCTCTCTACTTAGTCCAGCACGAAGTCTGGAGACCTCTGGCGGCAGCCTACCAAGAACAACCTGGACCGACCGGTGGTAC
CTCACCCCTTACCAGTCCGGCAGACAGTGTGGGTCCGCCGACACCAGACTAAGAACCTAGAACCTCGCTGGAAAAGGACCTTACAC
AGTCTGTCTGACCACCCCCACCGCCCTCAAAGTAGACGGCATCGCAGCTTGGATACACGCCGCCCACTGGCAGCAGAGTTTCC
CGACTGGAAAGCGGGCAGTGTAGCGCAACGCAATTAATGTGAGTTAGCTCCCAACTTTTAAAAAGAAAAGGGGGGATTGGGGGGTA
CAGTGCAGGGGAAAGAATAGTAGACATAATAGCAACAGACATACAACTAAAGAATTACAAAAACAATTTACAAAATTCAAAATTT
TTATCGATGCTTCCCGTCAACACCCCCCAACCCGCCCGGAGCTGAGAGTAATTCATACAAAAGGACTCGCCCTGCC
TTGGGGAATCCAGGGACCGTCTTAACTCCCACTAACGTAGAACCAGAGATCGCTGCGTTCCCGCCCCCTCACCCGCCCGCT
CTCGTCATCACTGAGGTGGAGAAGAGCATGCGTGAGGCTCCGGTGCCCGTCACTGGGCAGAGCGCACATCGCCACAGTCCCGA
GAAGTTGGGGGAGGGTTCGCAATTGAACCGGTGCTTAGAGAAGGTGGCGGGGTAACTGGGAAAGTGATGTCGTACTGG
CTCCGCTTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCCGTGAACGTTCTTTTTTCGCAACGGGTTTGCCG
CCAGAACACAGGTACACATATTGACCAAATCAGGGTAATTTTTGCATTTGTAATTTTTAAAAATGCTTTCTTTTAAATATACTT
TTTTGTTTATCTTATTTCTAATACTTTCCCTAATCTCTTTCTTTGAGGCAATAATGATACAATGTATCATGCCTCTTTGCACCA
TTCTAAAGAATAACAGTGATAATTTCTGGGTTAAGGCAATAGCGATTAGTTCTCGAGGATCCGACTGAAGTCGCTAGCTCGAC
TTTTGGAGAATATTTCTGCATATAAATATTTCTGCATATAAATTTGTAAGTCTGATGTAAGAGGTTTCATATTGCTAATAGCAGCTAC
AATCCAGCTACCATTCTGCTTTTTATTTTTATGGTTGGGATAAGGCTGGATTATTCTGAGTCCAAGCTAGGCCCTTTTGCTAATCAT
GTTTCATACCTCTTATCTTCTCCACAGCTCCTGGGCAACGTGCTGGTCTGTGTGCTGGCCATCACTTTGGCAAAGCACGTGAG
ATCTGAATTTCTGACACACCATGGTGAGCAAGGGCGAGGAGCTGTTACCCGGGTGGTGCCATCCTGGTCGAGCTGGACGGCGAC
GTAAACGGCCACAAGTTACAGCTGTCCGGCGAGGGCGAGGGCGATGCCACCTACGGCAAGCTGACCTGAAGTTCACTGACCA
CCGGCAAGCTGCCCCTGCCCACCCCTCGTGACCACCCCTGACCTACGGCGTGCAGTGTCTTACGCCCTACCCCGACCACAT
GAAGCAGCAGCTTCTTCAAGTCCGCCATGCCGAAGGCTACGTCCAGGAGCGCACCATCTTCTTCAAGGACGACGGCAACTAC
AAGCCCGCGCAGGTGAAGTTCGAGGGCAGCACCTTGGTGAACCGCATCGAGCTGAAGGGCATCGACTCAAGGAGGACGGCA
ACATCTGGGGCACAAGCTGGAGTACAACACTACAACAGCCACAACGCTTATATCATGCGCCGACAAGCAAGGCAAGGCAAGG
GAACTTCAAGATCCGCCACAACATCGAGGACGGCAGCGTGCAGCTCGCCGACCCTACCAGCAGAACACCCCATCGGCGACGGC
CCCGTGTCTGCTGCCGACAACCACTACCTGAGCACCAGTCCGCCCTGAGCAAAGACCCCAACGAGAAGCGCGATCACATGGTCC
TGCTGGAGTTCTGACCGCCCGGGATCACTCTCGGCATGGACGAGCTGTACAAGGCCACCATGACCGAGTACAAGCCACGGT
GCGCCTCGCCACCCGCGACGACGTCCCAGGGCCGTACGCACCCCTCGCCGCCCGCTTCGCCGACTACCCCGCCACGCGCCACACC
GTCGATCCGGACCGCCACATCGAGCGGGTACCGAGCTGCAAGAACTCTTCTCACGCGCGTCCGGCTCGACATCGGCAAGGTGT
GGGTTCGCGGACGACGGCGCCGCGGTGGCGGTCTGGACCACGCCGGAGAGCGTCAAGCGGGGGCGGTGTTTCGCCGAGATCGGCC
GCGATGGCCGAGTTGAGCGTTCCCGGTGGCCGCGCAGCAACAGATGGAAGGCTCCTGGCGCCGACCCGGCCCAAGGAGCCC
GCGTGGTTCTTGCCACCGTTCGGCGTCTCGCCGACCAACGCAAGGGTCTGGGACGCGCGTCTGCTCTCCCGGATGGAGG
CGGCCGAGCGCGCCGGGTGCCCGCTTCTTGAGACCTCCGCGCCCGCAACCTCCCTTCTACGAGCGGCTCGGCTTACCGT
CACCGCCGACGTGAGTGCCGAAGGACCGCGGACCTGGTGCATGACCCGCAAGCCCGGTGCTGAGCGGCCGACGACAGT
GTCGACGATAAAAATAAAGATTTTATTTAGTCTCCAGAAAAAGGGGGAAATGAAAGACCCACCTGTAGGTTTGGCAAGCTAGCT
TAAGTAACGCCATTTTGAAGGCATGAAAAATACATAACTGAGAATAGAGAAGTTGAGATCAAGGTCAGGAACAGATGGAACAG
CTGAATATGGGCCAAACAGGATATCTGTGGTAAGCAGTTTCTGCCCGGCTCAGGGCCAAGAACAGATGGAACAGCTGAATATGG
GCCAAACAGGATATCTGTGGTAAGCAGTTTCTGCCCGGCTCAGGGCCAAGAACAGATGGTCCCAGATGCGGTCCAGCCCTCAG
CAGTTTCTAGAGAACCATCAGATGTTTCCAGGGTGCCCAAGGACCTGAAATGACCTGTGCCTTATTTGAACCTAACCAATCAGT
TCGTTCTCGTTCTGTTGTCGCGCTTCTGCTCCCGAGCTCAATAAAAAGAGCCCAACCCCTCACTCGGGCGCCACTCTCC
GATTGACTGAGTCGCGCGGTACCCGTGTATCCAATAAACCTCTTTGACGTTGCATCCGACTGTGGTCTCGTGTCTTCTGGGA
GGGTCTCCTCTGAGTGATTGACTACCCGTGACGGGGTCTTTTCAATGCAGCATGTATCAAAAATAAATTTGGTTTTTTTTCTTA
AGTATTTACATTAATGACCATAGTTGCATTAATGAATCGCCAACGCGGGGAGAGGCGGTTTTCGTATTTGGCGCTCTTCCG
CTTCTCGTCACTGACTCGCTGCGTTCGGTCTGCGCTGCGCGAGCGGTATCAGTCACTCAAAGGCGGTAATACGGTTATC
CACAGAATCAGGGGATAACGCAGGAAAGAACATGTGAGCAAAAAGGCCAGCAAAAAGGCCAGGAACCGTAAAAAGGCCGCGTTGCTG
GCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAAACCCGACAGGACTAT
AAAGATAACAGGCGTTTCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCTGCCGCTTACCGGATACCTGTCCGCTT
TCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGTCAAGCTGTAGGTATCTCAGTTCGGTGTAGGTGTTTCGCTCCAAGCTGGGC
TGTGTGCACGAACCCCGTTACGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCCGTAAAGACAGACT
TATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCC
TAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGA AAAAGAGTTGGTAGTCTCT
TGATCCGGCAAAACAAACCACCGTGGTAGCGGTGGTTTTTTTTGTTTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAG
AAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGATTTTGGTTCATGAGATTATCAA
AAGGATCTTACCTAGATCTTTTTCGGGCCGGCCGAAATCAATCTAAAGTATATATGAGTAAACTTGGTCTGACAGTTACCAAT
GCTTAATCAGTGAGGCACCTATCTCAGCGATCTGTCTATTTTCGTTTATCCATAGTTGCCTGACTCCCCGTGCTGTAGATAACTAC
GATACGGGAGGGCTTACCATCTGGCCCCAGTGTGCAATGATACCGGAGACCCACGCTCACCGGCTCCAGATTATCAGCAATA
AACCAGCGAGCCGAAAGGGCCGAGCGAGAAGTGGTCTGCAACTTATCCGCTCCATCCAGTCTATTAATTTGTTGCCGGAAG
CTAGAGTAAGTAGTTGCCAGTTAATAAGTTTTCGCAACGTTGTTGCCATTGCTACAGGCATCGTGGTGTACGCTCGTCTGTTGG
TATGGCTTCAATCAGCTCCGGTTCCCAACGATCAAGGCGAGTTACATGATCCCCATGTTGTGCAAAAAAGCGGTTAGCTCCTTC
GGTCTCCGATCGTTGTCAGAAGTAAGTTGGCCGAGTGTATCACTCATGGTTATGGCAGCACTGCATAATTTCTTACTGTCA

TGCCATCCGTAAGATGCTTTTCTGTGACTGGTGAGTACTCAACCAAGTCATTCTGAGAATAGTGTATGCGGCGACCGAGTTGCTC
TTGCCCGGCGTCAATACGGGATAATACCGCGCCACATAGCAGAACTTTAAAAAGTGTCTCATCATTGGAAAAACGTTCTTCGGGGCGA
AAACTCTCAAGGATCTTACCGCTGTTGAGATCCAGTTCGATGTAACCCACTCGTGCACCCAACCTGATCTTCAGCATCTTTACTT
TCACCAGCGTTTCTGGGTGAGCAAAAAACAGGAAGGCAAAATGCCGCAAAAAAGGGAATAAGGGCGACACGGAAATGTTGAATACT
CATACTCTTCCTTTTTCAATATTATTGAAGCATTATCAGGGTTATTGTCTCATGAGCGGATACATATTTGAATGTATTTAGAAA
AATAAACAAATAGGGGTTCCGCGCACATTTT

License Information

pMXs vector is licensed from the University of Tokyo.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

This product is for RESEARCH USE ONLY; not for use in diagnostic procedures.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2012-2020: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.