

pMX- GFP Retroviral Vector

CATALOG NUMBER: RTV-050

STORAGE: -20°C

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

Background

Retroviruses are efficient tools for delivering heritable genes into the genome of dividing cells. Cell Biolabs' pMX-GFP retroviral vector is based on Moloney murine leukemia virus (MMLV). The vector provides the viral package signal, transcription and processing elements. 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*.

The vector contains the ampicillin-resistance gene, MMLV LTRs, package signal and GFP insert (Figure 1).

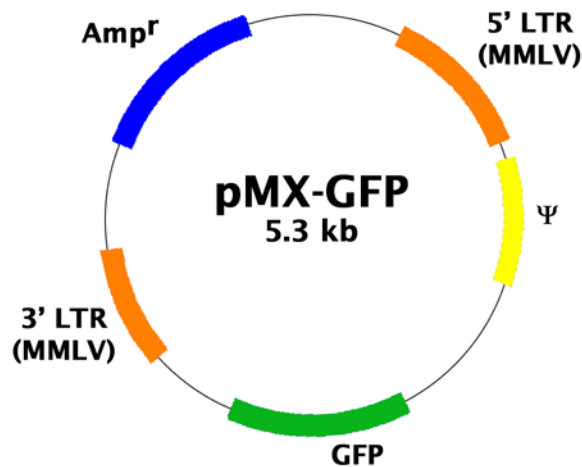


Figure 1. Schematic representation of pMX-GFP retroviral vector.

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.

References

1. Onishi, M., Kinoshita, S., Morikawa, Y., *et al.*, (1996) *Exp. Hematol.* **24**, 324-327.

Recent Production Citations

1. Verusingam, N.D., *et al.* (2017). Susceptibility of Human Oral Squamous Cell Carcinoma (OSCC) H103 and H376 cell lines to Retroviral OSKM mediated reprogramming. *PeerJ.* **5**:e3174. doi: 10.7717/peerj.3174. eCollection 2017.
2. Qu, Y. *et al.* (2016). NLRP3 recruitment by NLRC4 during Salmonella infection. *J Exp Med.* doi:10.1084/jem.20132234.
3. Malaver-Ortega, L. F. *et al.* (2015). Inducing pluripotency in cattle. *Methods Mol Biol.* doi:10.1007/978-1-4939-2848-4_6.
4. Parreno, J. *et al.* (2015). Efficient, low-cost nucleofection of passaged chondrocytes. *Cartilage.* doi:10.1177/1947603515609399.
5. Baird, A. E. G. *et al.* (2015). Derivation of canine induced pluripotent stem cells. *Reproduction in Domestic Animals.* DOI: 10.1111/rda.12562.
6. Gutiérrez-Fernández, A. *et al.* (2015). Loss of MT1-MMP causes cell senescence and nuclear defects which can be reversed by retinoic acid. *EMBO J.* doi:10.15252/embj.201490594.
7. Wattanapanitch, M. *et al.* (2014). Dual Small-Molecule Targeting of SMAD Signaling Stimulates Human Induced Pluripotent Stem Cells toward Neural Lineages *PLoS One.* **9**:e106952.
8. Gallaher, Z. R. *et al.* (2014). Neural proliferation in the dorsal root ganglia of the adult rat following capsaicin-induced neuronal death. *J Comp Neurol.* **522**:3295-3307.
9. Wahlestedt, M. *et al.* (2013). An epigenetic component of hematopoietic stem cell aging amenable to reprogramming into a young state. *Blood.* **121**:4257-4264.
10. Yi, L. *et al.* (2012). Multiple roles of p53-related pathways in somatic cell reprogramming and stem cell differentiation. *Cancer Res.* **72**: 5635-5645.

License Information

This product is licensed from Dr. Toshio Kitamura.

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

©2013-2017: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.