

pMXs-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' pMXs-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).

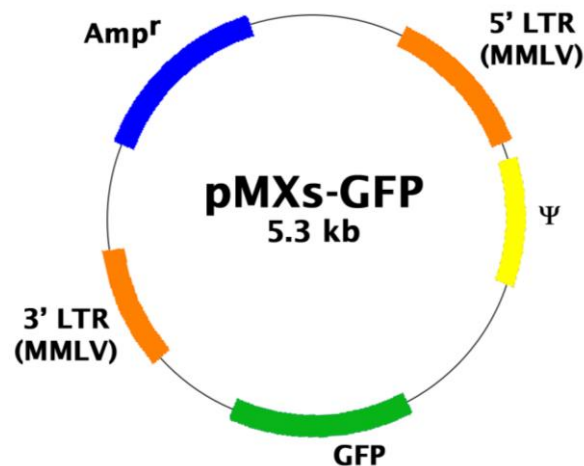


Figure 1. Schematic representation of pMXs-GFP retroviral vector.

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. Takeshita, M. *et al.* (2022). Potent SARS-CoV-2 neutralizing antibodies with therapeutic effects in two animal models. *iScience*. doi: 10.1016/j.isci.2022.105596.

2. Nagappan, A. et al. (2022). Caveolin-1-ACE2 axis modulates xenobiotic metabolism-linked chemoresistance in ovarian clear cell carcinoma. *Cell Biol Toxicol.* doi: 10.1007/s10565-022-09733-1.
3. Demirci, G. et al. (2021). (Bis)phosphonic Acid-Functionalized Poly(ethyleneimine)–Poly(amido amine)s for Selective In Vitro Transfection of Osteosarcoma Cells. *ACS Appl. Polym. Mater.* doi: 10.1021/acscapm.1c00297.
4. Nguyen, C. et al. (2021). Use of constitutive and inducible oncogene-containing iPSCs as surrogates for transgenic mice to study breast oncogenesis. *Stem Cell Res Ther.* **12**(1):301. doi: 10.1186/s13287-021-02285-x.
5. Kim, J. & Moon. Y. (2021). Mucosal ribosomal stress-induced PRDM1 promotes chemoresistance via stemness regulation. *Commun Biol.* **4**(1):543. doi: 10.1038/s42003-021-02078-1.
6. Petrova, S.C. et al. (2020). Regulation of breast cancer oncogenesis by the cell of origin's differentiation state. *Oncotarget.* **11**(43):3832-3848. doi: 10.18632/oncotarget.27783.
7. Kayagaki, N. et al. (2019). IRF2 transcriptionally induces GSDMD expression for pyroptosis. *Sci Signal.* **12**(582). pii: eaax4917. doi: 10.1126/scisignal.aax4917.
8. Terada, Y. et al. (2019). Human Pluripotent Stem Cell-Derived Tumor Model Uncovers the Embryonic Stem Cell Signature as a Key Driver in Atypical Teratoid/Rhabdoid Tumor. *Cell Rep.* **26**(10):2608-2621.e6. doi: 10.1016/j.celrep.2019.02.009.
9. Baird, A. et al. (2019). Biocompatible 3D printed thermoplastic scaffolds for osteoblast differentiation of equine iPS cells. *Tissue Eng Part C Methods.* doi: 10.1089/ten.TEC.2018.0343.
10. Thekkeparambil Chandrabose, S. et al. (2018). Amenable epigenetic traits of dental pulp stem cells underlie high capability of xeno-free episomal reprogramming. *Stem Cell Res Ther.* **9**(1):68. doi: 10.1186/s13287-018-0796-2.
11. Yakhkeshi, S. et al. (2018). In vitro improvement of quail primordial germ cell expansion through activation of TGF-beta signaling pathway. *J Cell Biochem.* **119**(6):4309-4319. doi: 10.1002/jcb.26618.
12. Takaya, T. et al. (2017). Autonomous xenogenic cell fusion of murine and chick skeletal muscle myoblasts. *Anim Sci J.* doi: 10.1111/asj.12884.
13. Kodaka, Y., Asakura, Y., and Asakura, A.(2017). Spin infection enables efficient gene delivery to muscle stem cells. *Biotechniques.* **63**(2):72-76. doi: 10.2144/000114576.
14. 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.
15. Qu, Y. et al. (2016). NLRP3 recruitment by NLRC4 during Salmonella infection. *J Exp Med.* doi:10.1084/jem.20132234.
16. Malaver-Ortega, L. F. et al. (2015). Inducing pluripotency in cattle. *Methods Mol Biol.* doi:10.1007/978-1-4939-2848-4_6.
17. Parreno, J. et al. (2015). Efficient, low-cost nucleofection of passaged chondrocytes. *Cartilage.* doi:10.1177/1947603515609399.
18. Baird, A. E. G. et al. (2015). Derivation of canine induced pluripotent stem cells. *Reproduction in Domestic Animals.* DOI: 10.1111/rda.12562.
19. 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.
20. 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.

21. 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.
22. Wahlestedt, M. et al. (2013). An epigenetic component of hematopoietic stem cell aging amenable to reprogramming into a young state. *Blood.* **121**:4257-4264.
23. 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.
5628 Copley Drive
San Diego, CA 92111
Worldwide: +1 858 271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2013-2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.