

pMYs-IRES-GFP Retroviral Vector

CATALOG NUMBER: RTV-021

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. Most retrovirus vectors including pBABE and pMXs are based on Moloney murine leukemia virus (MMLV). MMLV-based vectors usually are silenced in immature cells including embryonic carcinoma (EC) cells and embryonic stem (ES) cells, and possibly hematopoietic stem cells. Myeloproliferative sarcoma virus (MPSV) and PCC4-cell-passaged myeloproliferative sarcoma virus (PCMV) are mutants of MMLV and can stably express genes in immature cells including ES cells.

Cell Biolabs' pMYs-IRES-GFP retroviral vector (also known as pMYs-IG) includes hybrid LTRs containing elements from both MMLV and MPSV/PCMV, and it's capable of expressing genes in hematopoietic stem cells. 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*.

The vector contains the ampicillin-resistance gene, LTRs, package signal and MCS for cloning of your gene of interest (Figure 1).

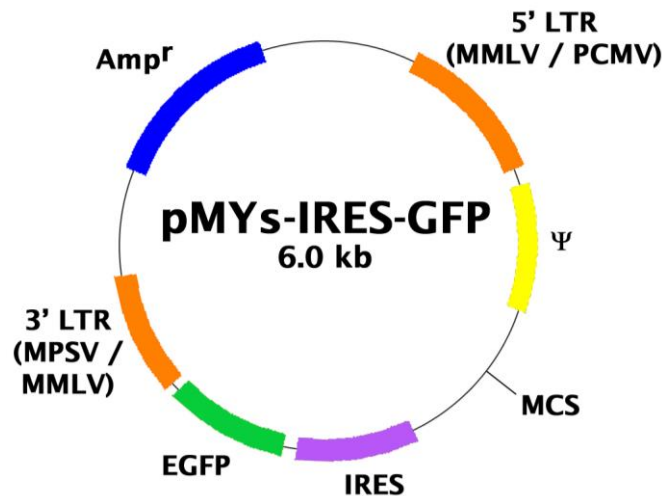


Figure 1. Schematic representation of pMYs-IRES-GFP retroviral vector.

MCS:

- Enzyme Sites: 5'-BamHI, EcoRI, XhoI, NotI, SnaBI-3'

- MCS Sequence:
TTAAGGATCCCAGTGTGGTGGTACGGGAATTCCTGCAGGCCTCGAGGGCCGGCGCGC
CGCGGCCGCTACGTAAATT---IRES---GFP---

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.

Reference

1. Kitamura T., *et al.*, (2003) *Exp. Hematol.* **31**, 1007-1014.

Recent Product Citations

1. Wang, G. *et al.* (2022). The RNA helicase DHX15 is a critical regulator of natural killer-cell homeostasis and functions. *Cell Mol Immunol.* doi: 10.1038/s41423-022-00852-7.
2. Gomaa, H.F. *et al.* (2022). Protective efficiency of Chelidonium majus extract against hepatotoxic and DNA changes induced by aflatoxin B1. *J. Appl. Pharm. Sci.* **12**(03): 140-149. doi: 10.7324/JAPS.2022.120315.
3. Uenaka, M. *et al.* (2022). Osteoblast-derived vesicles induce a switch from bone-formation to bone-resorption in vivo. *Nat Commun.* **13**(1):1066. doi: 10.1038/s41467-022-28673-2.
4. Xing, J. *et al.* (2021). Identification of poly(ADP-ribose) polymerase 9 (PARP9) as a noncanonical sensor for RNA virus in dendritic cells. *Nat Commun.* **12**(1):2681. doi: 10.1038/s41467-021-23003-4.
5. Murad, J.P. *et al.* (2021). Pre-conditioning Modifies the Tumor Microenvironment to Enhance Solid Tumor CAR T Cell Efficacy and Endogenous Protective Immunity. *Mol Ther.* doi: 10.1016/j.ymthe.2021.02.024.
6. Gaikwad, S.M. *et al.* (2020). A Small Molecule Stabilizer of the MYC G4-Quadruplex Induces Endoplasmic Reticulum Stress, Senescence and Pyroptosis in Multiple Myeloma. *Cancers (Basel).* **12**(10):E2952. doi: 10.3390/cancers12102952.
7. Muroy, S.E. *et al.* (2020). Phf15 - a novel transcriptional repressor regulating inflammation in a mouse microglial cell line. *Neuroimmunol Neuroinflammation.* doi: 10.20517/2347-8659.2020.16.
8. Zhang, S. *et al.* (2019). The transcription factor MZF1 differentially regulates murine Mtor promoter variants linked to tumor susceptibility. *J Biol Chem.* pii: jbc.RA119.009779. doi: 10.1074/jbc.RA119.009779.
9. 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.
10. Anandagoda, N. *et al.* (2019). microRNA-142-mediated repression of phosphodiesterase 3B critically regulates peripheral immune tolerance. *J Clin Invest.* **129**(3):1257-1271. doi: 10.1172/JCI124725.
11. Xia, Y. *et al.* (2019). The macrophage-specific V-ATPase subunit ATP6V0D2 restricts inflammasome activation and bacterial infection by facilitating autophagosome-lysosome fusion. *Autophagy.* 1-16. doi: 10.1080/15548627.2019.1569916.
12. Tanaka, M. *et al.* (2019). Development of a simple new flow cytometric antibody-dependent cellular cytotoxicity (ADCC) assay with excellent sensitivity. *J Immunol Methods.* **464**:74-86. doi: 10.1016/j.jim.2018.10.014.

13. Lin, Y. et al. (2018). Overexpression of Short Variant Form of New Kelch Family Protein Leads to Erythroid and Megakaryocyte Dysplasia by Targeting Megakaryocyte-Erythroid Progenitors. *DNA Cell Biol.* **37**(10):831-838. doi: 10.1089/dna.2018.4206.
14. Laurie, S.J. et al. (2018). 2B4 Mediates Inhibition of CD8+ T Cell Responses via Attenuation of Glycolysis and Cell Division. *J Immunol.* **201**(5):1536-1548. doi: 10.4049/jimmunol.1701240.
15. Zhang, M. et al. (2018). Transcription factor Hoxb5 reprograms B cells into functional T lymphocytes. *Nat Immunol.* **19**(3):279-290. doi: 10.1038/s41590-018-0046-x.
16. Weng, Q. et al. (2018). A protocol for generating induced T cells by reprogramming B cells in vivo. *Cell Regen (Lond).* **7**(1):7-15. doi: 10.1016/j.cr.2018.05.001.
17. Wang, T. et al. (2018). Trim27 confers myeloid hematopoiesis competitiveness by up-regulating myeloid master genes. *J Leukoc Biol.* **104**(4):799-809. doi: 10.1002/JLB.1A1217-480R.
18. Ye, B. et al. (2018). Klf4 glutamylation is required for cell reprogramming and early embryonic development in mice. *Nat Commun.* **9**(1):1261. doi: 10.1038/s41467-018-03008-2.
19. Humblin, E. et al. (2017). IRF8-dependent molecular complexes control the Th9 transcriptional program. *Nat Commun.* **8**(1):2085. doi: 10.1038/s41467-017-01070-w.
20. Liu, D. et al. (2016). Retrogenic ICOS expression increases differentiation of KLRG-1hiCD127loCD8+ T cells during Listeria infection and diminishes recall responses. *J Immunol.* doi:10.4049/jimmunol.1500218.
21. Xiao, X. et al. (2015). GITR subverts Foxp3+ Tregs to boost Th9 immunity through regulation of histone acetylation. *Nat Commun.* **6**:8266.
22. Chen, X. et al. (2015). OP9-Lhx2 stromal cells facilitate derivation of hematopoietic progenitors both in vitro and in vivo. *Stem Cell Res.* **15**:395-402.
23. Ogawara, Y. et al. (2015). IDH2 and NPM1 mutations cooperate to activate Hoxa9/Meis1 and hypoxia pathways in acute myeloid leukemia. *Cancer Res.* **75**:2005-2016.
24. Amin, S. et al. (2015). Hoxa2 selectively enhances meis binding to change a branchial arch ground state. *Dev Cell.* **9**:265-277.
25. Yang, D. et al. (2015). Enforced expression of hoxa5 in hematopoietic stem cells leads to aberrant erythropoiesis in vivo. *Cell Cycle.* doi:10.4161/15384101.2014.992191.
26. Zvezdova, E. et al. (2014). In vivo functional mapping of the conserved protein domains within murine Themis1. *Immunol Cell Biol.* **92**:721-728.
27. Ye, B. et al. (2014). Cytosolic carboxypeptidase CCP6 is required for megakaryopoiesis by modulating Mad2 polyglutamylation. *J Exp Med.* **11**:2439-2454.
28. Baba, T. et al. (2012). Novel process of intrathymic tumor-immune tolerance through CCR2-mediated recruitment of Sirpα+ dendritic cells: a murine model. *PLoS One.* **7**(7):e41154. doi: 10.1371/journal.pone.0041154.

License Information

This product 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' 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

©2012-2024: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.