

pMXs-IRES-GFP Retroviral Vector

CATALOG NUMBER: RTV-013

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-IRES-GFP retroviral vector (also known as pMXs-IG) 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*.

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

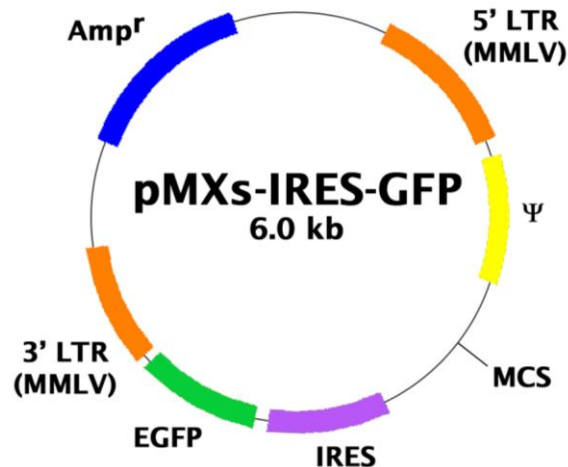


Figure 1. Schematic representation of pMXs-IRES-GFP retroviral vector.

MCS:

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

TTAATTAAGGATCCCCAGTGTGGTGGTACGGGAATTCCTGCAGGCCTCGAGGGCCGGC
GCGCCGCGCCGCTACGTAAATT---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. Kusano, S. et al. (2023). Contributions of the N-terminal flanking residues of an antigenic peptide from the Japanese cedar pollen allergen Cry j 1 to the T-cell activation by HLA-DP5. *Int Immunol.* doi: 10.1093/intimm/dxad024.
2. Wu, A. et al. (2023). Structural basis for the allosteric modulation of rhodopsin by nanobody binding to its extracellular domain. *Nat Commun.* **14**(1):5209. doi: 10.1038/s41467-023-40911-9.
3. Okada, M. et al. (2023). Detection of mutant antigen-specific T cell receptor against multiple myeloma for T cell engineering. *Mol Ther Methods Clin Dev.* doi: 10.1016/j.omtm.2023.05.014.
4. Zwicklhuber, J. et al. (2023). A Novel Fluorescence-Based Screen of Gene Editing Molecules for Junctional Epidermolysis Bullosa. *Int J Mol Sci.* **24**(6):5197. doi: 10.3390/ijms24065197.
5. Kobayashi, E. et al. (2023). Gene modified NK cell line as a powerful tool for evaluation of cloned TCRs for TCR-T cell therapy. *Cell Immunol.* doi: 10.1016/j.cellimm.2022.104656.
6. Tsukumo, S.I. et al. (2022). AFF3, a susceptibility factor for autoimmune diseases, is a molecular facilitator of immunoglobulin class switch recombination. *Sci Adv.* **8**(34):eabq0008. doi: 10.1126/sciadv.abq0008.
7. Ogawa, H. et al. (2022). Nectin-2 Acts as a Viral Entry Mediated Molecule That Binds to Human Herpesvirus 6B Glycoprotein B. *Viruses.* **14**(1):160. doi: 10.3390/v14010160.
8. Hwang, S.J. et al. (2022). miR-125a-5p attenuates macrophage-mediated vascular dysfunction by targeting Ninjurin1. *Cell Death Differ.* doi: 10.1038/s41418-021-00911-y.
9. Yamaguchi, S. et al. (2021). TCR function analysis using a novel system reveals the multiple unconventional tumor-reactive T cells in human breast cancer-infiltrating lymphocytes. *Eur J Immunol.* doi: 10.1002/eji.202049070.
10. Bhat, S. et al. (2021). ZNF471 modulates EMT and functions as methylation regulated tumor suppressor with diagnostic and prognostic significance in cervical cancer. *Cell Biol Toxicol.* doi: 10.1007/s10565-021-09582-4.
11. Deng, Z. et al. (2020). Def6 regulates endogenous type-I interferon responses in osteoblasts and suppresses osteogenesis. *Elife.* **9**:e59659. doi: 10.7554/eLife.59659.
12. Mou, Z. et al. (2020). Identification of a Protective Leishmania Antigen Dihydrolipoyl Dehydrogenase and Its Responding CD4+ T Cells at Clonal Level. *J Immunol.* doi: 10.4049/jimmunol.2000338.
13. Bourdely, P. et al. (2020). Transcriptional and Functional Analysis of CD1c+ Human Dendritic Cells Identifies a CD163+ Subset Priming CD8+CD103+ T Cells. *Immunity.* S1074-7613(20)30232-6. doi: 10.1016/j.immuni.2020.06.002.
14. Cai, Z. et al. (2020). ROR agonist hampers the proliferation and survival of post-activated CD8+ T cells through the downregulation of cholesterol synthesis-related genes. *Immunol Cell Biol.* doi: 10.1111/imcb.12406.
15. Yamakawa, T. et al. (2020). Novel gene Merlot inhibits differentiation and promotes apoptosis of osteoclasts. *Bone.* doi: 10.1016/j.bone.2020.115494.
16. Mashima, H. et al. (2020). The role of calcium-binding protein S100g (CalbindinD-9K) and annexin A10 in acute pancreatitis. *Biochem Biophys Res Commun.* pii: S0006-291X(20)30653-7. doi: 10.1016/j.bbrc.2020.03.155.
17. Fujiwara, K. et al. (2020). Impact of scFv structure in chimeric antigen receptor on receptor expression efficiency and antigen recognition properties. *Biochem Biophys Res Commun.* pii: S0006-291X(20)30556-8. doi: 10.1016/j.bbrc.2020.03.071.
18. Mutoh, T. et al. (2020). Identification of two major autoantigens negatively regulating endothelial activation in Takayasu arteritis. *Nat Commun.* **11**(1):1253. doi: 10.1038/s41467-020-15088-0.

19. Fujinami, H. et al. (2020). CLP1 acts as the main RNA kinase in mice. *Biochem Biophys Res Commun.* pii: S0006-291X(20)30334-X. doi: 10.1016/j.bbrc.2020.02.066.
20. Cella, M. et al. (2019). Subsets of ILC3-ILC1-like cells generate a diversity spectrum of innate lymphoid cells in human mucosal tissues. *Nat Immunol.* doi: 10.1038/s41590-019-0425-y.
21. Barrow, A.D. et al. (2018). Natural Killer Cells Control Tumor Growth by Sensing a Growth Factor. *Cell.* **172**(3):534-548.e19. doi: 10.1016/j.cell.2017.11.037.
22. Methot, S.P. et al. (2018). A licensing step links AID to transcription elongation for mutagenesis in B cells. *Nat Commun.* **9**(1):1248. doi: 10.1038/s41467-018-03387-6.
23. Avbelj, M. et al. (2018). The role of N-terminal segment and membrane association in MyD88-mediated signaling. *Biochem Biophys Res Commun.* **495**(1):878-883. doi: 10.1016/j.bbrc.2017.11.099.
24. Jung, H.J. (2017). The Ubiquitin-like with PHD and Ring Finger Domains 1 (UHRF1)/DNA Methyltransferase 1 (DNMT1) Axis is a Primary Regulator of Cell Senescence. *J Biol Chem.* **292**(9):3729-3739. doi: 10.1074/jbc.M116.750539.
25. Mueller, A.A. et al. (2016). Intronic polyadenylation of PDGF α in resident stem cells attenuates muscle fibrosis. *Nature* **540**:276-279.
26. Baek, J.M. et al. (2016). Nicotinamide phosphoribosyltransferase inhibits receptor activator of nuclear factor- κ B ligand-induced osteoclast differentiation in vitro. *Mol. Med. Rep.* doi:10.3892/mmr.2016.6069.
27. Chiu, Y-H. et al. (2016). Dendritic cell-specific transmembrane protein (DC-STAMP) regulates osteoclast differentiation via the Ca²⁺/NFATc1 axis. *J. Cellular Physiol.* 10.1002/jcp.25638.
28. Kimura, T. et al. (2016). Polarization of M2 macrophages requires Lamtor1 that integrates cytokine and amino acid signals. *Nature Comm.* doi:10.1038/ncomms13130.
29. Werner, S. et al. (2015). Iroquois homeobox 2 suppresses cellular motility and chemokine expression in breast cancer cells. *BMC Cancer.* **15**:896.
30. Ravikumar, P. et al. (2015). Nanoparticle facilitated inhalational delivery of erythropoietin receptor cDNA protects against hyperoxic lung injury. *Nanomedicine.* doi:10.1016/j.nano.2015.10.004.

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'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

©2008-2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.