

## pMXs- IRES- GFP Retroviral Vector

**CATALOG NUMBER:** RTV-013

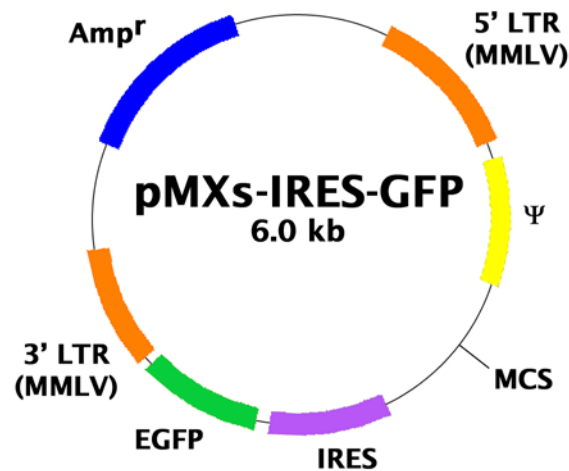
**STORAGE:** -20°C

**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:

TTAATTAAGGATCCCAGTGTGGTGGTACGGGAATTCCTGCAGGCCTCGAGGGCCGGC  
 GCGCCGCGGCCGCTACGTAAATT---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.

## References

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

## Recent Product Citations

1. 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.
2. Mueller, A.A. et al. (2016). Intronic polyadenylation of PDGF $\alpha$  in resident stem cells attenuates muscle fibrosis. *Nature* **540**:276-279.
3. Baek, J.M. et al. (2016). Nicotinamide phosphoribosyltransferase inhibits receptor activator of nuclear factor-kB ligand-induced osteoclast differentiation in vitro. *Mol. Med. Rep.* doi:10.3892/mmr.2016.6069.
4. Chiu, Y-H. et al. (2016). Dendritic cell-specific transmembrane protein (DC-STAMP) regulates osteoclast differentiation via the Ca<sup>2+</sup>/NFATc1 axis. *J. Cellular Physiol.* 10.1002/jcp.25638.
5. Kimura, T. et al. (2016). Polarization of M2 macrophages requires Lamtor1 that integrates cytokine and amino acid signals. *Nature Comm.* doi:10.1038/ncomms13130.
6. Werner, S. et al. (2015). Iroquois homeobox 2 suppresses cellular motility and chemokine expression in breast cancer cells. *BMC Cancer.* **15**:896.
7. 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.
8. Amagai, Y. et al. (2015). A point mutation in the extracellular domain of KIT promotes tumorigenesis of mast cells via ligand-independent auto-dimerization. *Sci Rep.* **5**:9775.
9. Sakata-Yanagimoto, M. & Chiba, S. (2015). Notch2 signaling in mast cell development and distribution in the intestine. *Methods Mol Biol.* **1220**:79-89.
10. Shapiro, I.M. et al. (2014). Merlin deficiency predicts FAK inhibitor sensitivity: a synthetic lethal relationship. *Sci Transl Med.* **237**:237ra68.
11. Zhang, Q. et al. (2013). TNF- $\alpha$  impairs differentiation and function of TGF- $\beta$ -induced Treg cells in autoimmune diseases through Akt and Smad3 signaling pathway. *J Mol Cell Biol.* 10.1093/jmcb/mjs063.
12. Koso, H. et al. (2012). PNAS Plus: Transposon mutagenesis identifies genes that transform neural stem cells into glioma-initiating cells. *PNAS.* 10.1073/pnas.1215899109.
13. Cui, G. et al. (2009). Berberine differentially modulates the activities of ERK, p38 MAPK, and JNK to suppress Th17 and Th1 T cell differentiation in Type 1 diabetic mice. *J. Biol. Chem.* **284**:28420-28429.
14. Guan, Y. et al. (2010). Human TLRs 10 and 1 share common mechanisms of innate immune sensing but not signaling. *J. Immunol.* **184**:5094-5103.

## 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.  
7758 Arjons Drive  
San Diego, CA 92126  
Worldwide: +1 858-271-6500  
USA Toll-Free: 1-888-CBL-0505  
E-mail: [tech@cellbiolabs.com](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2008-2017: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.