

## pMXs-IRES-Blasticidin Retroviral Vector

**CATALOG NUMBER:** RTV-016

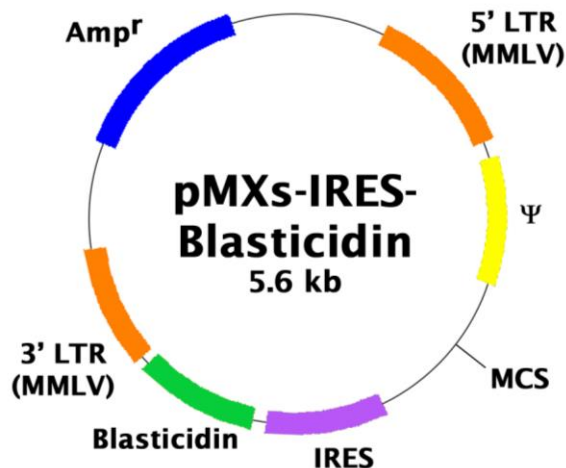
**STORAGE:** -20°C

**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-Blasticidin retroviral vector (also known as pMXs-IB) 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-Blasticidin retroviral vector.

MCS:

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

### 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

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

## Recent Product Citations

1. Mizuike, A. et al. (2023). The C10orf76-PI4KB axis orchestrates CERT-mediated ceramide trafficking to the distal Golgi. *J Cell Biol.* **222**(7):e202111069. doi: 10.1083/jcb.202111069.
2. Laqtom, N.N. et al. (2022). CLN3 is required for the clearance of glycerophosphodiester from lysosomes. *Nature.* **609**(7929):1005-1011. doi: 10.1038/s41586-022-05221-y.
3. Krzyzanowska, A.K. et al. (2021). Activation of nuclear factor-kappa B by TNF promotes nucleus pulposus mineralization through inhibition of ANKH and ENPP1. *Sci Rep.* **11**(1):8271. doi: 10.1038/s41598-021-87665-2.
4. Zhu, X.G. et al. (2020). Functional Genomics In Vivo Reveal Metabolic Dependencies of Pancreatic Cancer Cells. *Cell Metab.* doi: 10.1016/j.cmet.2020.10.017.
5. Sviderskiy, V.O. et al. (2020). Hyperactive CDK2 Activity in Basal-like Breast Cancer Imposes a Genome Integrity Liability that Can Be Exploited by Targeting DNA Polymerase  $\epsilon$ . *Mol Cell.* doi: 10.1016/j.molcel.2020.10.016.
6. Watanabe, M. et al. (2020). A substrate-trapping strategy to find E3 ubiquitin ligase substrates identifies Parkin and TRIM28 targets. *Commun Biol.* **3**(1):592. doi: 10.1038/s42003-020-01328-y.
7. Wimmer, M. et al. (2020). A cancer stem cell-like phenotype is associated with miR-10b expression in aggressive squamous cell carcinomas. *Cell Commun Signal.* **18**(1):61. doi: 10.1186/s12964-020-00550-9.
8. Williams, R.T. et al. (2020). ZBTB1 Regulates Asparagine Synthesis and Leukemia Cell Response to L-Asparaginase. *Cell Metab.* **31**(4):852-861.e6. doi: 10.1016/j.cmet.2020.03.008.
9. Tamura, I. et al. (2020). Wilms tumor 1 regulates lipid accumulation in human endometrial stromal cells during decidualization. *J Biol Chem.* pii: jbc.RA120.012841. doi: 10.1074/jbc.RA120.012841.
10. Wang, C. et al. (2019). PARP1 hinders histone H2B occupancy at the NFATc1 promoter to restrain osteoclast differentiation. *J Bone Miner Res.* doi: 10.1002/jbmr.3927.
11. Musiani, D. et al. (2019). Proteomics profiling of arginine methylation defines PRMT5 substrate specificity. *Sci Signal.* **12**(575). pii: eaat8388. doi: 10.1126/scisignal.aat8388.
12. Peng, W. et al. (2019). Functional analysis of the secondary HIV-1 capsid binding site in the host protein cyclophilin A. *Retrovirology.* **16**(1):10. doi: 10.1186/s12977-019-0471-4.
13. Zhu, X.G. et al. (2019). CHP1 Regulates Compartmentalized Glycerolipid Synthesis by Activating GPAT4. *Mol Cell.* **74**(1):45-58.e7. doi: 10.1016/j.molcel.2019.01.037.
14. Bunse, L. et al. (2018). Suppression of antitumor T cell immunity by the oncometabolite (R)-2-hydroxyglutarate. *Nat Med.* **24**(8):1192-1203. doi: 10.1038/s41591-018-0095-6.
15. Dovey, C.M. et al. (2018). MLKL Requires the Inositol Phosphate Code to Execute Necroptosis. *Mol Cell.* **70**(5):936-948.e7. doi: 10.1016/j.molcel.2018.05.010.
16. Ganier, O. et al. (2018). Structural centrosome aberrations sensitize polarized epithelia to basal cell extrusion. *Open Biol.* **8**(6). pii: 180044. doi: 10.1098/rsob.180044.
17. Qi, C. et al. (2018). Anti-mitotic chemotherapeutics promote apoptosis through TL1A-activated death receptor 3 in cancer cells. *Cell Res.* **28**(5):544-555. doi: 10.1038/s41422-018-0018-6.
18. Take, K., et al. (2017). CDK5 Regulatory Subunit-Associated Protein 1-like 1 Negatively Regulates Adipocyte Differentiation through Activation of Wnt Signaling Pathway. *Sci Rep.* **7**(1):7326. doi: 10.1038/s41598-017-06469-5.

19. Law, M. E. et al. (2016). CUB domain-containing protein 1 and the epidermal growth factor receptor cooperate to induce cell detachment. *Breast Cancer Res.* **18**:80.
20. Saito, A. et al. (2016). Roles of capsid-interacting host factors in multimodal inhibition of HIV-1 by PF74. *Jf Virol.* doi:10.1128/JVI.03116-15.
21. Peking, P. et al. (2016). A gene gun-mediated nonviral RNA trans-splicing strategy for Col7a1 repair. *Mol Ther Nucleic Acids.* doi:10.1038/mtna.2016.3.
22. Sotillo, E. et al. (2015). Convergence of acquired mutations and alternative splicing of CD19 enables resistance to CART-19 immunotherapy. *Cancer Discov.* **5**:1282-1295.
23. Greil, C. et al. (2015). The role of APC/CCdh1 in replication stress and origin of genomic instability. *Oncogene.* doi:10.1038/onc.2015.367.
24. Choi, Y. A. et al. (2014). Bobby Sox homology regulates odontoblast differentiation of human dental pulp stem cells/progenitors. *Cell Commun Signal.* **12**:35.
25. Papatheodorou, P. et al. (2011). Lipolysis-stimulated lipoprotein receptor (LSR) is the host receptor for the binary toxin clostridium difficile transferase (CDT). *PNAS* **108**: 16422-16427.
26. Sugatani, T. et al. (2011). A microRNA expression signature of osteoclastogenesis. *Blood* **117**:3648-3657.

### **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](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2008-2024: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.