

pMXs-IRES-Neo Retroviral Vector

CATALOG NUMBER: RTV-015

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-Neo retroviral vector (also known as pMXs-IN) 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).

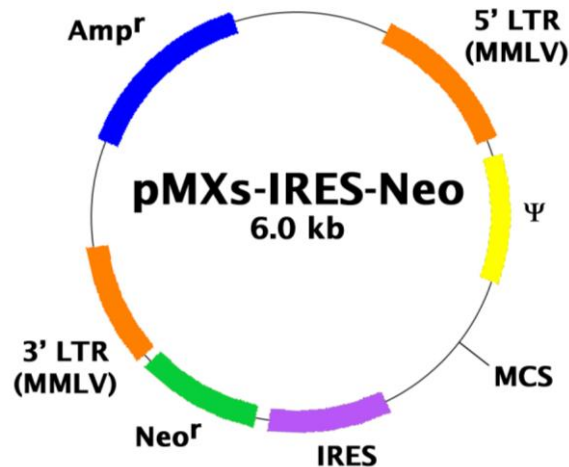


Figure 1. Schematic representation of pMXs-IRES-Neo retroviral vector.

MCS:

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

```
TTAATTAAGGATCCCAGTGTGGTGGTACGGGAATTCCTGCAGGCCTCGAGGGCCGGC
GCGCCGCGGCCGCTACGTAAATT---IRES---neo---
```

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. da Silva Almeida, A. *et al.* (2025). Building a potent TREM2 agonistic, biparatopic, common light chain antibody. *MAbs*. **17**(1):2546554. doi: 10.1080/19420862.2025.2546554.
2. Hassenrück, F. *et al.* (2023). Functional impact and molecular binding modes of drugs that target the PI3K isoform p110 δ . *Commun Biol.* **6**(1):603. doi: 10.1038/s42003-023-04921-z.
3. Goto, A. *et al.* (2022). Compartmentalization of casein kinase 1 γ CSNK1G controls the intracellular trafficking of ceramide. *iScience*. doi: 10.1016/j.isci.2022.104624.
4. Maemura, T. *et al.* (2021). Antibody-Dependent Enhancement of SARS-CoV-2 Infection Is Mediated by the IgG Receptors Fc γ RIIA and Fc γ RIIIA but Does Not Contribute to Aberrant Cytokine Production by Macrophages. *mBio*. doi: 10.1128/mBio.01987-21.
5. 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.
6. Kuroda, M. *et al.* (2020). Identification of interferon-stimulated genes that attenuate Ebola virus infection. *Nat Commun.* **11**(1):2953. doi: 10.1038/s41467-020-16768-7.
7. Mashima, H. *et al.* (2018). The role of G α q/G α 11 signaling in intestinal epithelial cells. *Biochem Biophys Res.* **13**:93-98. doi: 10.1016/j.bbrep.2018.01.003.
8. Koso, H. *et al.* (2014). Identification of FoxR2 as an oncogene in medulloblastoma. *Cancer Res.* **74**:2351-2361.

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.
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

©2012-2025: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.