

## pBABE-Puro Retroviral Vector

**CATALOG NUMBER:** RTV-001-PURO

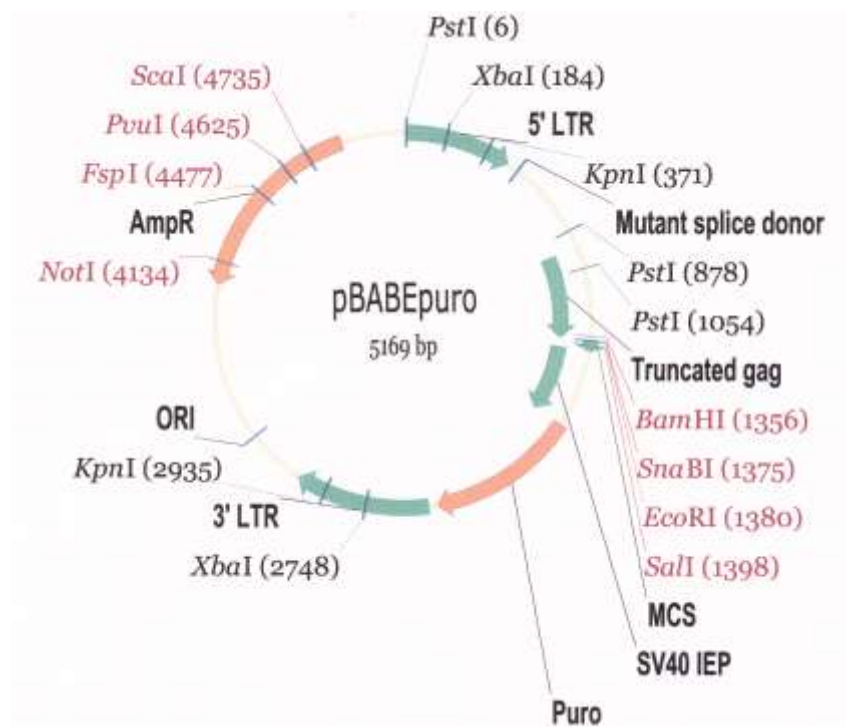
**STORAGE:** -20°C

**QUANTITY AND CONCENTRATION:** 10 µg at 0.5 µg/µL in TE

### Background

Retroviruses are efficient tools for delivering heritable genes into the genome of dividing cells. Cell Biolabs’ retrovirus vector is based on the pBABE vector system, which is derived from Moloney murine leukemia virus (MMLV). The vector provides the viral package signal, transcription and processing elements, and 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 are used to express silencing RNAs (siRNA) to decrease the expression of target genes both *in vitro* and *in vivo*.

The vector contains the bacterial origin of replication, ampicillin-resistance gene, and puromycin-resistance gene for the growth of infected mammalian cells to select stable cell lines (Figure 1).



**Figure 1.** pBABE-Puro Retroviral Vector Map

## **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. Morgenstern, J. P. and H Land. (1990) *Nuc. Acid Res.* 18, 3587-3596.
2. Coffin, J. M. and H. E. Varmus, *Retroviruses*, Cold Spring Harbor Press, NY.
3. Schuck S, Manninen A, Honsho M, Fullekrug J and Simons K. (2004) *Proc Natl Acad Sci U S A.* 101, 4912-4917.

## **Recent Product Citations**

1. Zhou, H.C. et al. (2021). Dual and opposing roles of the androgen receptor in VETC-dependent and invasion-dependent metastasis of hepatocellular carcinoma. *J Hepatol.* doi: 10.1016/j.jhep.2021.04.053.
2. Handa, O. et al. (2021). PDZK1 induces resistance to apoptosis in esophageal adenocarcinoma cells. *Esophagus.* doi: 10.1007/s10388-021-00819-z.
3. Yang, C. et al. (2021). Nuclear IGF1R interacts with NuMA and regulates 53BP1-dependent DNA double-strand break repair in colorectal cancer. *Oncol Rep.* **46**(2):168. doi: 10.3892/or.2021.8119.
4. Keenan, S.N. et al. (2020). Perilipin 5 S155 phosphorylation by PKA is required for the control of hepatic lipid metabolism and glycemic control. *J Lipid Res.* doi: 10.1194/jlr.RA120001126.
5. Wu, K.Z.L. et al. (2019). Pathogenic FAM83G palmoplantar keratoderma mutations inhibit the PAWS1:CK1 $\alpha$  association and attenuate Wnt signalling. *Wellcome Open Res.* **4**:133. doi: 10.12688/wellcomeopenres.15403.1.
6. Wenta, T. et al. (2019). HtrA4 Protease Promotes Chemotherapeutic-Dependent Cancer Cell Death. *Cells.* **8**(10):1112. doi: 10.3390/cells8101112.
7. Hatanaka, T. et al. (2019). The neural ELAVL protein HuB enhances endogenous proto-oncogene activation. *Biochem Biophys Res Commun.* pii: S0006-291X(19)31443-3. doi: 10.1016/j.bbrc.2019.07.089.
8. Wenta, T. et al. (2019). The HtrA3 protease promotes drug-induced death of lung cancer cells by cleavage of the X-linked inhibitor of apoptosis protein (XIAP). *FEBS J.* doi: 10.1111/febs.14977.
9. Katamune, C. et al. (2019). Mutation of the gene encoding the circadian clock component PERIOD2 in oncogenic cells confers chemoresistance by up-regulating the Aldh3a1 gene. *J Biol Chem.* **294**(2):547-558. doi: 10.1074/jbc.RA118.004942.
10. Sang, Y. et al. (2018). Snail promotes metastasis of nasopharyngeal carcinoma partly by down-regulating TEL2. *Cancer Commun (Lond).* **38**(1):58. doi: 10.1186/s40880-018-0328-6.
11. Hu, H. et al. (2018). Long non-coding RNA XLOC\_000647 suppresses progression of pancreatic cancer and decreases epithelial-mesenchymal transition-induced cell invasion by down-regulating NLRP3. *Mol Cancer.* **17**(1):18. doi: 10.1186/s12943-018-0761-9.
12. Fulcher, L.J. et al. (2017). Targeting endogenous proteins for degradation through the affinity-directed protein missile system. *Open Biol.* doi:10.1098/rsob.170066.
13. Dong, G. et al. (2017). Crk adaptor proteins regulate CD3 $\zeta$  chain phosphorylation and TCR/CD3 down-modulation in activated T cells. *Cell Signal.* **36**:117-126.
14. Ji, X. et al. (2017). Permissive, in vitro replication of hepatitis B virus genotype E. *J Virol Methods.* **243**:20-24. doi: 10.1016/j.jviromet.2017.01.016.

15. Lin, Y. et al. (2017). SUMO-modified Insulin-Like Growth Factor 1 Receptor (IGF-1R) Increases Cell Cycle Progression and Cell Proliferation. *J Cell Physiol*. doi: 10.1002/jcp.25818.
16. Jiang, J. et al. (2016). Overexpression of microRNA-21 strengthens stem cell-like characteristics in a hepatocellular carcinoma cell line. *World J. Surg. Onc.* **14**: 278.
17. Fulcher, L.J. et al. (2016). An affinity-directed protein missile system for targeted proteolysis. *Open Biol.* 10.1098/rsob.160255.
18. Na, H. N. et al. (2016). E4orf1 enhances glucose uptake independent of proximal insulin signaling. *PLoS One.* **11**:e0161275.
19. Kohda, A. et al. (2016). The nuclear protein IκBζ forms a transcriptionally active complex with nuclear factor-κB (NF-κB) p50 and Lcn2 promoter via the N-and C-terminal ankyrin repeat motifs. *J Biol Chem*. doi:10.1074/jbc.M116.719302.
20. Mathur, P. S. et al. (2016). Kinase-dependent and-independent roles for PTK6 in colon cancer. *Mol Cancer Res*. doi:10.1158/1541-7786.MCR-15-0450.
21. Challita-Eid, P. M. et al. (2016). Enfortumab vedotin antibody–drug conjugate targeting nectin-4 is a highly potent therapeutic agent in multiple preclinical cancer models. *Cancer Res.* **76**:3003-3013.
22. Verhoeven, R. J. et al. (2016). NF-κB signaling regulates expression of Epstein-Barr virus BART microRNAs and long noncoding RNAs in nasopharyngeal carcinoma. *J Virol*. doi:10.1128/JVI.00613-16.
23. Endo, S. et al. (2016). TEX101, a glycoprotein essential for sperm fertility, is required for stable expression of Ly6k on testicular germ cells. *Sci Rep*. doi:10.1038/srep23616.
24. Hegde, V. et al. (2015). An Adenovirus-derived protein: A novel candidate for anti-diabetic drug development. *Biochimie*. doi:10.1016/j.biochi.2015.12.002.
25. Rai, R. et al. (2015). Arginyltransferase suppresses cell tumorigenic potential and inversely correlates with metastases in human cancers. *Oncogene*. doi:10.1038/onc.2015.473.
26. Matveeva, E. A. et al. (2015). Vimentin is involved in regulation of mitochondrial motility and membrane potential by Rac1. *Biol Open*. doi: 10.1242/bio.011874.
27. Chakroun, I. et al. (2015). Genome-wide association between Six4, MyoD and the histone demethylase Utx during myogenesis. *FASEB J*. doi:10.1096/fj.15-277053.
28. Hansen, M. T. et al. (2015). A link between inflammation and metastasis: serum amyloid A1 and A3 induce metastasis, and are targets of metastasis-inducing S100A4. *Oncogene*. **34**:424-435.
29. Wei, M. et al. (2015). The over-expression of aquaporin-1 alters erythroid gene expression in human erythroleukemia K562 cells. *Tumour Biol*. **36**:1-12.
30. Jiao, H. K. et al. (2015). Prognostic significance of Cbx4 expression and its beneficial effect for transarterial chemoembolization in hepatocellular carcinoma. *Cell Death Disease*. **6**:e1689.

### **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)

©2007-2022: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.