

## Platinum-A Retroviral Packaging Cell Line, Amphotropic

**CATALOG NUMBER:** RV-102

**STORAGE:** Liquid nitrogen

*Note: For best results begin culture of cells immediately upon receipt. If this is not possible, store at -80°C until first culture. Store subsequent cultured cells long term in liquid nitrogen.*

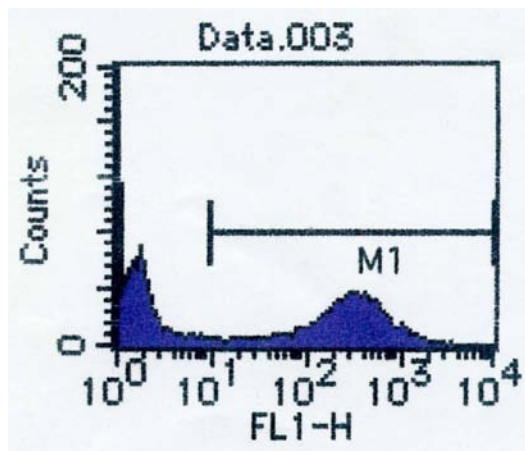
**QUANTITY & CONCENTRATION:** 1.0 mL,  $>3 \times 10^6$  cells/mL in DMEM, 20% FCS and 10% DMSO

### **Background**

Retroviruses are efficient tools for delivering heritable genes into the genome of dividing cells. However, conventional NIH-3T3 based retroviral packaging cell lines have limited stability and produce low viral yields, mainly due to poor expression level of the retroviral structure proteins (gag, pol, env) in the packaging cells.

The Platinum-A (Plat-A) Cell Line, a potent retrovirus packaging cell line based on the 293T cell line, was generated using novel packaging constructs with an EF1 $\alpha$  promoter to ensure longer stability and high-yield retroviral structure protein expression (gag, pol, amphotropic env). Plat-A cells can be kept in good condition in for at least 4 months in the presence of drug selection, and can produce retroviruses with an average titer of  $1 \times 10^6$  infectious units/mL by transient transfection. In addition, replication competent retroviruses (RCR) are virtually nonexistent because only coding sequences of viral structural genes are used, avoiding any unnecessary retroviral sequences.

The Plat-A cell line is designed for rapid, transient production of high-titer, amphotropic retrovirus.



**Figure 1.** TF-1 cells were infected with GFP retrovirus supernatant produced in Plat-A cells after transfection with pMX-GFP.

## Medium

1. Culture Medium: DMEM, 10% fetal calf serum (FCS), 1  $\mu\text{g}/\text{mL}$  puromycin, 10  $\mu\text{g}/\text{mL}$  blasticidin, penicillin and streptomycin
2. Freeze Medium: 70% DMEM, 20% FBS, 10% DMSO

## Methods

### **I. Establishing Plat-A Cultures from Frozen Cells**

1. After quickly thawing the cells in a 37°C water bath, immediately transfer the thawed cell suspension into a 15 mL tube containing 10 mL of culture medium.
2. Centrifuge the tube for 5 min at 1300 to 1500 rpm.
3. Discard the supernatant and break the cell pellet by finger tapping.
4. Add a few drops of culture medium with gentle shaking and finger tap the tube a few times.
5. Add 2 mL of culture medium to the tube and gently pipet the cell suspension up and down twice.
6. Transfer the cell suspension to a 10 cm culture dish (Falcon® #3003 works well) containing 8 mL of culture medium.
7. Swirl the culture plate well to mix the cells, then incubate the cells for three days before expansion.

#### ***Important Notes:***

- ***Don't change the culture medium during the first three days. It is normal to see some cells floating after the first 24 hours.***
- ***Don't culture cells to complete confluency. Split cells 4X to 6X every two to three days when the culture reaches 70-90% confluency.***

### **II. Splitting the Cells**

*Note: Avoid forming bubbles as much as possible during this procedure.*

1. Wash cells once with PBS.
2. Add 4 mL of 0.05% Trypsin/0.5 mM EDTA solution to a 10 cm dish and incubate at 37°C for 3-5 min.
3. Remove the cells from the dish surface by tapping the rim of the culture dish.
4. Transfer 10 mL of the culture medium to a 50 mL tube.
5. Using the same pipette with some residual culture medium, wash the dish surface gently three times in 4 mL of the Trypsin/EDTA solution.
6. Gently pipette the cell suspension up and down 7 times and transfer the cell suspension into the 50 mL tube containing 10 mL medium from step 4.
7. Centrifuge the cells for 5 min at 1300-1500 rpm.
8. Discard the supernatant and break the cell pellet by finger tapping.
9. Add a few drops of culture medium with gentle shaking and finger tap the tube a few times.
10. Add 5 mL of culture medium and gently pipet the cell suspension up and down twice.
11. Add 15 mL of culture medium, then count and seed the cells. Typically  $10^7$  cells can be harvested from one 10 cm culture dish.

## Transfection

1. Seed  $2 \times 10^6$  cells in a 60 mm culture dish without antibiotics including puromycin and blasticidin one day before transfection.
2. After 16 to 24 hours, start transfection when the culture becomes 70-80% confluent.  
*Note: We suggest transfecting cells with FuGENE® Transfection Reagent (Roche Applied Science) or Lipofectamine™ Plus (Invitrogen). For example, 3 µg retroviral expression plasmid is mixed with 9 µL FuGENE® Transfection Reagent according to the manufacturer's recommendation. The mixed DNA- FuGENE® complex is added by dropwise into the culture media.*
3. Harvest retroviral supernatant 48 hours after transfection.

## References

1. Morita, S., Kojim, T., and Kitamura, T. (2000) *Gene Therapy* 7: 1063-1066.

## Recent Product Citations

1. Grasso S, et al. (2017). The scaffold protein p140Cap limits ERBB2-mediated breast cancer progression interfering with Rac GTPase-controlled circuitries. *Nat Commun.* **8**:14797. doi: 10.1038/ncomms14797.
2. Hu, T. et al. (2016). Integrin  $\alpha 6$ /Akt/Erk signaling is essential for human breast cancer resistance to radiotherapy. *Sci Rep.* doi:10.1038/srep33376.
3. Patel, P. L. et al. (2016). Derepression of hTERT gene expression promotes escape from oncogene-induced cellular senescence. *Proc Natl Acad Sci U S A.* doi:10.1073/pnas.1602379113.
4. Ha, Y. J. et al. (2016). SOCS1 suppresses IL-1 $\beta$ -induced C/EBP $\beta$  expression via transcriptional regulation in human chondrocytes. *Exp Mol Med.* doi:10.1038/emmm.2016.47.
5. Huang, C. et al. (2016). SUMOylated ORC2 recruits a histone demethylase to regulate centromeric histone modification and genomic stability. *Cell Rep.* doi:10.1016/j.celrep.2016.02.091.
6. Koseki, J. et al. (2016). A trans-omics mathematical analysis reveals novel functions of the ornithine metabolic pathway in cancer stem cells. *Sci Rep.* **6**:20726.
7. Naemura, M. et al. (2015). ANRIL regulates the proliferation of human colorectal cancer cells in both two-and three-dimensional culture. *Mol Cell Biochem.* doi:10.1007/s11010-015-2618-5.
8. Friedrich Ben-Nun, I. et al. (2015). Generation of induced pluripotent stem cells from mammalian endangered species. *Methods Mol Biol.* doi:10.1007/978-1-4939-2848-4\_10 101-109.
9. Massouridès, E. et al. (2015). Dp412e: a novel human embryonic dystrophin isoform induced by BMP4 in early differentiated cells. *Skelet Muscle.* **5**:40.
10. Chinyenetere, F. et al. (2015). Mice null for the deubiquitinase USP18 spontaneously develop leiomyosarcomas. *BMC Cancer.* **15**:886.
11. Guan, H. et al. (2015). FOXO1 inhibits osteosarcoma oncogenesis via Wnt/ $\beta$ -catenin pathway suppression. *Oncogenesis.* doi:10.1038/oncsis.2015.25.
12. Chen, Q. et al. (2015). Chitosan prevents adhesion during rabbit flexor tendon repair via the sirtuin 1 signaling pathway. *Mol Med Rep.* **12**:4598-4603.
13. Chien, Y. et al. (2015). Synergistic effects of carboxymethyl-hexanoyl chitosan, cationic polyurethane-short branch PEI in miR122 gene delivery: Accelerated differentiation of iPSCs into mature hepatocyte-like cells and improved stem cell therapy in a hepatic failure model. *Acta Biomater.* **13**:228-244.

14. Fuerstenau-Sharp, M. et al. (2015). Generation of highly purified human cardiomyocytes from peripheral blood mononuclear cell-derived induced pluripotent stem cells. *PLoS One*. **10**:e0126596.
15. Kwon, H. J. et al. (2015). NK cell function triggered by multiple activating receptors is negatively regulated by glycogen synthase kinase-3 $\beta$ . *Cell Signal*. doi:10.1016/j.cellsig.2015.05.012.
16. Lee, J. H. et al. (2015). AKT phosphorylates H3-threonine 45 to facilitate termination of gene transcription in response to DNA damage. *Nucleic Acids Res*. doi: 10.1093/nar/gkv176.
17. Han, F. et al. (2015). Human induced pluripotent stem cell-derived neurons improve motor asymmetry in a 6-hydroxydopamine-induced rat model of Parkinson's disease. *Cytotherapy*. doi: 10.1016/j.jcyt.2015.02.001.
18. Balboni, A. L. et al. (2015). p53 and DeltaNp63 $\alpha$  co-regulate the transcriptional and cellular response to TGF $\beta$  and BMP Signals. *Mol Cancer Res*. doi: 10.1158/1541-7786.MCR-14-0152-T.
19. Ji, Y. et al. (2015). miR-155 augments CD8+ t-cell antitumor activity in lymphoreplete hosts by enhancing responsiveness to homeostatic  $\gamma$ C Cytokines. *Proc Natl Acad Sci U S A*. **112**:476-481.
20. DeCastro, A. J. et al. (2015).  $\Delta$ NP63 $\alpha$  transcriptionally activates chemokine receptor 4 (CXCR4) expression to regulate breast cancer stem cell activity and chemotaxis. *Mol Cancer Ther*. **14**:225-235.
21. Hu, R. et al. (2015). NF- $\kappa$ B signaling is required for XBP1 (unspliced and spliced)-mediated effects on antiestrogen responsiveness and cell fate decisions in breast cancer. *Mol Cell Biol*. **35**:379-390.
22. Robin, J. D. et al. (2015). Isolation and immortalization of patient-derived cell lines from muscle biopsy for disease modeling. *J Vis Exp*. **95**:e52307-e52307.
23. Lui, V. W. et al. (2014). Frequent mutation of receptor protein tyrosine phosphatases provides a mechanism for STAT3 hyperactivation in head and neck cancer. *Proc Natl Acad Sci U S A*. **111**:1114-1119.
24. Qian, H. et al. (2014). PCH-2 regulates *Caenorhabditis elegans* lifespan. *Aging*. **7**.
25. Wong, C. et al. (2014). The bromodomain and extra-terminal inhibitor CPI203 enhances the antiproliferative effects of rapamycin on human neuroendocrine tumors. *Cell Death Dis*. **5**:e1450.
26. Goh, P. A. & Verma, P. J. (2014). Generation of induced pluripotent stem cells from mouse adipose tissue. *Methods Mol Biol*. **1194**:253-270.
27. Kondoh, T. et al. (2014). CD7 promotes extramedullary involvement of the B-cell acute lymphoblastic leukemia line Tanoue by enhancing integrin  $\beta$ 2-dependent cell adhesiveness. *Int J Oncol*. **45**:1073-1081.
28. Wattanapanitch, M. et al. (2014). Dual Small-Molecule Targeting of SMAD Signaling Stimulates Human Induced Pluripotent Stem Cells toward Neural Lineages *PLoS One*. **9**:e106952.
29. Hayashi, K. et al. (2014). Visualization and characterization of cancer stem-like cells in cervical cancer. *Int J Oncol*. **45**:2468-2474.
30. Hamabe, A. et al. (2014). Role of pyruvate kinase M2 in transcriptional regulation leading to epithelial-mesenchymal transition. *Proc Natl Acad Sci U S A*. **111**:15526-15531.
31. Krivega, I. et al. (2014). Role of LDB1 in the transition from chromatin looping to transcription activation. *Genes Dev*. **28**:1279-1290.

## **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.

## **License Information**

This licensed product is intended for ACADEMIC, GOVERNMENT AND NON-PROFIT RESEARCH USE ONLY; not for use in diagnostic or therapeutic procedures. The product may not be transferred, sold, or otherwise provided to another laboratory except by an authorized distributor of Cell Biolabs, Inc.

Use of this product by Biotechnology and Pharmaceutical companies requires a license for all fields of use including research. Please contact:

Director of Business Development  
Cell Biolabs, Inc.  
[busdev@cellbiolabs.com](mailto:busdev@cellbiolabs.com)

## **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.