

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 X 10⁶ cells/mL in DMEM, 20% FBS 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 α 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 for at least 4 months in the presence of drug selection, and can produce retroviruses with an average titer of 1 x 10⁶ 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.

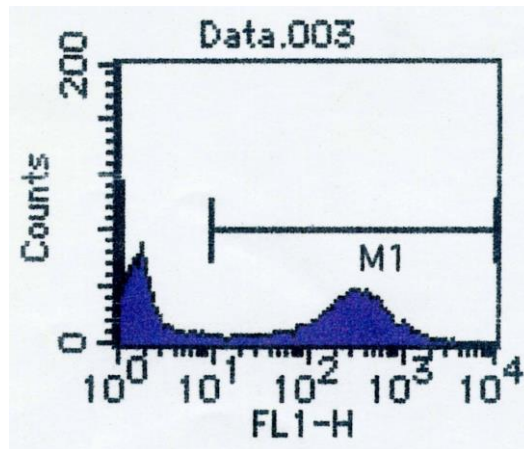


Figure 1. TF-1 cells were infected with GFP retrovirus supernatant produced in Plat-A cells after transfection with pMX-GFP.

Medium

1. Culture Medium: DMEM, 10% fetal bovine serum (FBS), 1 μ g/mL puromycin, 10 μ g/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 2×10^6 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.

Reference

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

Recent Product Citations

1. Rojas-Restrepo, J. et al. (2023). Functional Relevance of CTLA4 Variants: an Upgraded Approach to Assess CTLA4-Dependent Transendocytosis by Flow Cytometry. *J Clin Immunol*. doi: 10.1007/s10875-023-01582-9.
2. Sako, H. et al. (2023). microRNAs slow translating ribosomes to prevent protein misfolding in eukaryotes. *EMBO J*. doi: 10.15252/embj.2022112469.
3. Cannell, I.G. et al. (2023). FOXC2 promotes vasculogenic mimicry and resistance to anti-angiogenic therapy. *Cell Rep*. **42**(8):112791. doi: 10.1016/j.celrep.2023.112791.
4. Harada, Y. et al. (2023). Metabolic clogging of mannose triggers dNTP loss and genomic instability in human cancer cells. *Elife*. **12**:e83870. doi: 10.7554/eLife.83870.
5. Torcal Garcia, G. et al. (2023). Carm1-arginine methylation of the transcription factor C/EBP α regulates transdifferentiation velocity. *Elife*. doi: 10.7554/eLife.83951.
6. Lei, Y. et al. (2023). Cooperative sensing of mitochondrial DNA by ZBP1 and cGAS promotes cardiotoxicity. *Cell*. doi: 10.1016/j.cell.2023.05.039.
7. Cho, M.K. et al. (2023). Frankincense ameliorates endometriosis via inducing apoptosis and reducing adhesion. *Integr Med Res*. **12**(2):100947. doi: 10.1016/j.imr.2023.100947.
8. Wu, S.J. et al. (2023). Immunotherapeutic potential of blinatumomab-secreting γ 9 δ 2 T Cells. *Transl Oncol*. doi: 10.1016/j.tranon.2023.101650.
9. Miura, K. et al. (2023). Chorioallantoic membrane assay revealed the role of TIPARP (2,3,7,8-tetrachlorodibenzo-p-dioxin-inducible poly (ADP-ribose) polymerase) in lung adenocarcinoma-induced angiogenesis. *Cancer Cell Int*. **23**(1):34. doi: 10.1186/s12935-023-02870-5.
10. Zenke, S. et al. (2022). Differential trafficking of ligands trogocytosed via CD28 versus CTLA4 promotes collective cellular control of co-stimulation. *Nat Commun*. **13**(1):6459. doi: 10.1038/s41467-022-34156-1.
11. Ueda, Y. et al. (2022). Mechanistic insights into cancer drug resistance through optogenetic PI3K signaling hyperactivation. *Cell Chem Biol*. doi: 10.1016/j.chembiol.2022.10.002.
12. Choi, S. et al. (2022). Enhanced tumor targeting and timely viral release of mesenchymal stem cells/oncolytic virus complex due to GRP78 and inducible E1B55K expressions greatly increase the antitumor effect of systemic treatment. *Mol Ther Oncolytics*. doi: 10.1016/j.omto.2022.09.004.
13. Masuta, Y. et al. (2022). Assessment of Fc γ receptor-dependent binding of influenza hemagglutinin vaccine-induced antibodies in a non-human primate model. *iScience*. **25**(10):105085. doi: 10.1016/j.isci.2022.105085.
14. Sakakibara, M. et al. (2022). Bitter taste receptor T2R38 is expressed on skin-infiltrating lymphocytes and regulates lymphocyte migration. *Sci Rep*. **12**(1):11790. doi: 10.1038/s41598-022-15999-6.
15. Seo, S.H. et al. (2022). PTEN/AKT signaling pathway related to hTERT downregulation and telomere shortening induced in Toxoplasma GRA16-expressing colorectal cancer cells. *Biomed Pharmacother*. doi: 10.1016/j.biopha.2022.113366.
16. Shih, H.T. et al. (2022). DNMT3b protects centromere integrity by restricting R-loop-mediated DNA damage. *Cell Death Dis*. **13**(6):546. doi: 10.1038/s41419-022-04989-1.

17. Kapadia, B.B. et al. (2022). PARK2 regulates eIF4B-driven lymphomagenesis. *Mol Cancer Res.* doi: 10.1158/1541-7786.MCR-21-0729.
18. Kim, H.J. et al. (2021). SLAC2B-dependent microtubule acetylation regulates extracellular matrix-mediated intracellular TM4SF5 traffic to the plasma membranes. *FASEB J.* **35**(3):e21369. doi: 10.1096/fj.202002138RR.
19. Yarmarkovich, M. et al. (2021). Cross-HLA targeting of intracellular oncoproteins with peptide-centric CARs. *Nature.* **599**(7885):477-484. doi: 10.1038/s41586-021-04061-6.
20. Gao, R. et al. (2021). YAP/TAZ and ATF4 drive resistance to Sorafenib in hepatocellular carcinoma by preventing ferroptosis. *EMBO Mol Med.* doi: 10.15252/emmm.202114351.
21. Bai, T. et al. (2021). Establishment of human induced trophoblast stem-like cells from term villous cytotrophoblasts. *Stem Cell Res.* **56**:102507. doi: 10.1016/j.scr.2021.102507.
22. Mellis, I.A. et al. (2021). Responsiveness to perturbations is a hallmark of transcription factors that maintain cell identity in vitro. *Cell Syst.* **12**(9):885-899.e8. doi: 10.1016/j.cels.2021.07.003.
23. Lesch, S. et al. (2021). T cells armed with C-X-C chemokine receptor type 6 enhance adoptive cell therapy for pancreatic tumours. *Nat Biomed Eng.* doi: 10.1038/s41551-021-00737-6.
24. 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.
25. Kim, J. & Moon, Y. (2021). Mucosal ribosomal stress-induced PRDM1 promotes chemoresistance via stemness regulation. *Commun Biol.* **4**(1):543. doi: 10.1038/s42003-021-02078-1.
26. Hyun, Y.M. et al. (2020). Endogenous DEL-1 restrains melanoma lung metastasis by limiting myeloid cell-associated lung inflammation. *Sci Adv.* **6**(45):eabc4882. doi: 10.1126/sciadv.abc4882.
27. Petrova, S.C. et al. (2020). Regulation of breast cancer oncogenesis by the cell of origin's differentiation state. *Oncotarget.* **11**(43):3832-3848. doi: 10.18632/oncotarget.27783.
28. Nitta, S. et al. (2020). Three-dimensional spheroid culture of canine hepatocyte-like cells derived from bone marrow mesenchymal stem cells. *Regen Ther.* doi: 10.1016/j.reth.2020.09.002.
29. Luanpitpong, S. et al. (2020). A novel TRPM7/O-GlcNAc axis mediates tumour cell motility and metastasis by stabilising c-Myc and caveolin-1 in lung carcinoma. *Br J Cancer.* doi: 10.1038/s41416-020-0991-7.
30. Putri, D.S. et al. (2020). Cytomegalovirus infection elicits a conserved chemokine response from human and Guinea pig amnion cells. *Virology.* doi: 10.1016/j.virol.2020.06.005.

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

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

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