

## A549/GFP Cell Line

**CATALOG NUMBER:** AKR-209

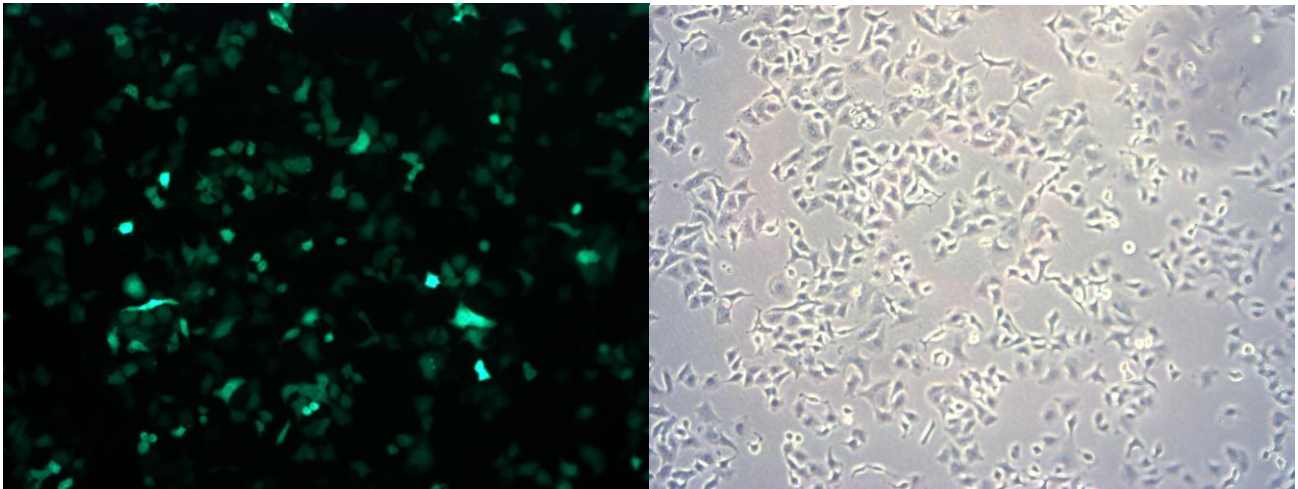
**STORAGE:** Liquid nitrogen

*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 mL,  $1 \times 10^6$  cells/mL in 70% DMEM, 20% FBS, 10% DMSO

### **Background**

The A549 cell line was first developed in 1972 through the removal and culturing of cancerous lung tissue in the explanted tumor of 58-year-old male. A549 cells are human alveolar basal epithelial cells. They are squamous in nature and responsible for the diffusion of substances across the alveoli of lungs. In vitro they grow adherently as a monolayer, and in vivo they induce tumors in athymic mice. A549 cells could synthesize lecithin with a high percentage of desaturated fatty acids utilizing the cytidine diphosphocholine pathway. Our A549/GFP cell line stably expresses GFP; the gene was introduced using lentivirus.



**Figure 1.** A549/GFP Cell Line. Left: GFP Fluorescence; Right: Phase Contrast.

### **Quality Control**

This cryovial contains at least  $1.0 \times 10^6$  A549/GFP cells as determined by morphology, trypan-blue dye exclusion, and viable cell count. The A549/GFP cells are tested free of microbial contamination.

### **Medium**

1. Culture Medium: D-MEM (high glucose), 10% fetal bovine serum (FBS), 0.1 mM MEM Non-Essential Amino Acids (NEAA), 2 mM L-glutamine, 1% Pen-Strep.
2. Freeze Medium: 70% DMEM, 20% FBS, 10% DMSO.

## **Methods**

### **Establishing A549/GFP Cultures from Frozen Cells**

1. Place 10 mL of complete DMEM or McCoy's 5A growth medium in a 50-mL conical tube. Thaw the frozen cryovial of cells within 1–2 minutes by gentle agitation in a 37°C water bath. Decontaminate the cryovial by wiping the surface of the vial with 70% (v/v) ethanol.
2. Transfer the thawed cell suspension to the conical tube containing 10 ml of growth medium.
3. Collect the cells by centrifugation at 1000 rpm for 5 minutes at room temperature. Remove the growth medium by aspiration.
4. Resuspend the cells in the conical tube in 15 mL of fresh growth medium by gently pipetting up and down.
5. Transfer the 15 mL of cell suspension to a T-75 tissue culture flask. Place the cells in a 37°C incubator at 5% CO<sub>2</sub>.
6. Monitor cell density daily. Cells should be passaged when the culture reaches 95% confluence.

### **Recent Product Citations**

1. Mora, D.S.O. et al. (2023). An optimized live bacterial delivery vehicle safely and efficaciously delivers bacterially transcribed therapeutic nucleic acids. *Eng Life Sci.* **23**(3): e2200037. doi: 10.1002/elsc.202200037.
2. Mendieta, I. et al. (2021). Ultrastructural changes associated to the neuroendocrine transdifferentiation of the lung adenocarcinoma cell line A549. *Acta Histochem.* **123**(8):151797. doi: 10.1016/j.acthis.2021.151797.
3. Wen, Y. et al. (2020). A supramolecular platform for controlling and optimizing molecular architectures of siRNA targeted delivery vehicles. *Sci Adv.* **6**(31):eabc2148. doi: 10.1126/sciadv.abc2148.
4. Ni, B.S. et al. (2019). Plug-and-Play In Vitro Metastasis System toward Recapitulating the Metastatic Cascade. *Sci Rep.* **9**(1):18110. doi: 10.1038/s41598-019-54711-z.
5. Obeid, M.A. et al. (2017). Formulation of nonionic surfactant vesicles (NISV) prepared by microfluidics for therapeutic delivery of siRNA into cancer cells. *Mol. Pharm.* **14**(7):2450-2458.
6. Kumar, A. et al. (2017). Influenza virus exploits tunneling nanotubes for cell-to-cell spread. *Scientific Reports.* **7**: 40360. doi:10.1038/srep40360
7. Shopsowitz, K. E. et al. (2015). Periodic-shRNA molecules are capable of gene silencing, cytotoxicity and innate immune activation in cancer cells. *Nucleic Acids Res.* doi:10.1093/nar/gkv1488.
8. Almosnid, N. M. et al. (2015). In vitro antitumor effects of two novel oligostilbenes, cis-and trans-suffruticosol D, isolated from *Paeonia suffruticosa* seeds. *Int J Oncol.* doi:10.3892/ijo.2015.3269.
9. Zhu, L. et al. (2014). Matrix metalloproteinase 2-sensitive multifunctional polymeric micelles for tumor-specific co-delivery of siRNA and hydrophobic drugs. *Biomaterials.* **35**:4213-4222.

### **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)

©2010-2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.