

293AAV Cell Line

CATALOG NUMBER: AAV-100

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 x 10⁶ cells/mL in 90% complete medium, 10% DMSO

Background

Adeno-associated virus (AAV) belongs to the family of Parvoviridae, a group of viruses among the smallest of single-stranded and non-enveloped DNA viruses. There are nine different AAV serotypes reported to date. AAV can infect both dividing and non-dividing cells and can be maintained in the human host cell, creating the potential for long-term gene transfer. Recombinant AAV-2 is the most common serotype used in gene delivery, and can be produced at high titers with a helper virus or Cell Biolabs' AAV Helper-Free System in 293 cells.

The 293AAV Cell Line is a permanent line established from primary embryonic human kidney transformed with human adenovirus type 5 DNA. The genes encoded by the E1 region of adenovirus (E1a and E1b) are expressed in these cells and participate in transactivation of viral promoters, allowing these cells to produce high levels of protein.

293AAV is derived from the parental 293 cell line, through cloning and multiple rounds of testing, 293AAV is specifically selected for a high level of AAV production in a helper-free system. It offers several advantages over the regular 293 cells:

- Larger cell surface area resulting higher transfection and better yield of AAV.
- Flattened morphology.
- Firm attachment to culture plate, ideal for large scale culture and AAV production.

Quality Control

This cryovial contains at least 1.0 × 10⁶ 293AAV cells as determined by morphology, trypan-blue dye exclusion, and viable cell count. The 293AAV 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 (optional)
2. Freeze Medium: 90% complete medium, 10% DMSO

Methods

Establishing 293AAV Cultures from Frozen Cells

1. Place 10 mL of complete DMEM 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% CO₂.
6. Monitor cell density daily. Cells should be passaged when the culture reaches 95% confluence.

Recent Product Citations

1. Yun, T. et al. (2021). Inhibitor of DNA binding 2 (Id2) mediates microtubule polymerization in the brain by regulating α K40 acetylation of α -tubulin. *Cell Death Discov.* **7**(1):257. doi: 10.1038/s41420-021-00652-4.
2. Chao, G. et al. (2021). Measurement of Large Serine Integrase Enzymatic Characteristics in HEK293 Cells Reveals Variability and Influence on Downstream Reporter Expression. *FEBS J.* doi: 10.1111/febs.16037.
3. Sveidahl Johansen, O. et al. (2021). Lipolysis drives expression of the constitutively active receptor GPR3 to induce adipose thermogenesis. *Cell.* doi: 10.1016/j.cell.2021.04.037.
4. Hu, J. et al. (2021). Co-opting regulation bypass repair as a gene correction strategy for monogenic diseases. *Mol Ther.* doi: 10.1016/j.ymthe.2021.04.017.
5. Motori, E. et al. (2020). Larsson, Neuronal metabolic rewiring promotes resilience to neurodegeneration caused by mitochondrial dysfunction. *Sci. Adv.* **6**(35). doi: 10.1126/sciadv.aba8271.
6. Göbel, J. et al. (2020). Mitochondria-Endoplasmic Reticulum Contacts in Reactive Astrocytes Promote Vascular Remodeling. *Cell Metab.* pii: S1550-4131(20)30120-0. doi: 10.1016/j.cmet.2020.03.005.
7. Lin, Y.H. et al. (2020). Accumbal D2R-medium spiny neurons regulate aversive behaviors through PKA-Rap1 pathway. *Neurochem Int.* doi: 10.1016/j.neuint.2020.104935.
8. Siegrist, C.M. et al. (2020). CRISPR/Cas9 as an antiviral against Orthopoxviruses using an AAV vector. *Sci Rep.* **10**(1):19307. doi: 10.1038/s41598-020-76449-9.
9. Sun, W. et al. (2020). snRNA-seq reveals a subpopulation of adipocytes that regulates thermogenesis. *Nature.* **587**(7832):98-102. doi: 10.1038/s41586-020-2856-x.
10. Bougioukli, S. et al. (2020). Limited potential of AAV-mediated gene therapy in transducing human mesenchymal stem cells for bone repair applications. *Gene Ther.* doi: 10.1038/s41434-020-0182-4.
11. Chen, S.H. et al. (2020). A Simple, Two-Step, Small-Scale Purification of Recombinant Adeno-Associated Viruses. *J Virol Methods.* doi: 10.1016/j.jviromet.2020.113863.
12. Wu, Z. et al. (2020). Gene therapy conversion of striatal astrocytes into GABAergic neurons in mouse models of Huntington's disease. *Nat Commun.* **11**(1):1105. doi: 10.1038/s41467-020-14855-3.

13. Miyawaki, T. et al. (2020). Visualization and molecular characterization of whole-brain vascular networks with capillary resolution. *Nat Commun.* **11**(1):1104. doi: 10.1038/s41467-020-14786-z.
14. Funahashi, Y. et al. (2019). Phosphorylation of Npas4 by MAPK Regulates Reward-Related Gene Expression and Behaviors. *Cell Rep.* **29**(10):3235-3252.e9. doi: 10.1016/j.celrep.2019.10.116.
15. Nakano, R. et al. (2019). Auxin-mediated rapid degradation of target proteins in hippocampal neurons. *Neuroreport.* **30**(13):908-913. doi: 10.1097/WNR.0000000000001299.
16. Kuzumaki, N. et al. (2019). Cell-specific overexpression of COMT in dopaminergic neurons of Parkinson's disease. *Brain.* doi:10.1093/brain/awz084.
17. Ao, H. et al. (2019). Egr1 mediates retinal vascular dysfunction in diabetes mellitus via promoting p53 transcription. *J Cell Mol Med.* doi: 10.1111/jcmm.14225.
18. Nakamura, M. et al. (2019). Glycogen Synthase Kinase-3 α Promotes Fatty Acid Uptake and Lipotoxic Cardiomyopathy. *Cell Metab.* pii: S1550-4131(19)30005-1. doi: 10.1016/j.cmet.2019.01.005.
19. Iwasawa, C. et al. (2019). Regional and temporal regulation and role of somatostatin receptor subtypes in the mouse brain following systemic kainate-induced acute seizures. *Neurosci Res.* pii: S0168-0102(18)30670-9. doi: 10.1016/j.neures.2019.01.004.
20. Hashimoto, H. et al. (2016). Study on AAV-mediated gene therapy for diabetes in humanized liver mouse to predict efficacy in humans. *Biochem Biophys Res Commun.* doi:10.1016/j.bbrc.2016.08.104.
21. Lee, S. E. et al. (2016). nArgBP2 regulates excitatory synapse formation by controlling dendritic spine morphology. *Proc Natl Acad Sci U S A.* doi:10.1073/pnas.1600944113.
22. Lutz, D. et al. (2014). Myelin basic protein cleaves cell adhesion molecule L1 and promotes neurogenesis and cell survival. *J Biol Chem.* **289**:13503-13518.

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
www.cellbiolabs.com

©2009-2021: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.

