

293 LTV Cell Line

CATALOG NUMBER: LTV-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, $>1 \times 10^6$ cells/mL in 90% complete medium, 10% DMSO

Background

The 293LTV 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.

293LTV also stably expresses the SV40 large T antigen and Neomycin resistance gene, through cloning and multiple rounds of testing for viral yield, 293LTV is specifically selected for high level of lentiviral production. It offers several advantages over the regular 293T cells:

- High lentiviral yield
- Firm attachment to culture plate and fast growing
- Ideal as a host when making lentivirus by cotransfection

Quality Control

This cryovial contains at least 1.0×10^6 293LTV cells as determined by morphology, trypan-blue dye exclusion, and viable cell count. The 293LTV 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

I. Establishing 293LTV 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. Wade, M. et al. (2021). In vivo generation of collagen specific Tregs with AAV8 suppresses autoimmune responses and arthritis in DBA1 mice through IL10 production. *Sci Rep.* **11**(1):18204. doi: 10.1038/s41598-021-97739-w.
2. Smith, C.M. et al. (2021). Phosphatidylserine binding directly regulates TIM-3 function. *Biochem J.* **478**(17):3331-3349. doi: 10.1042/BCJ20210425.
3. Kiss, E. et al. (2021). Artesunate restores the levels of inhibitory synapse proteins and reduces amyloid- β and C-terminal fragments (CTFs) of the amyloid precursor protein in an AD-mouse model. *Mol Cell Neurosci.* doi: 10.1016/j.mcn.2021.103624.
4. Kiss, E. et al. (2021). Artemisinin-treatment in pre-symptomatic APP-PS1 mice increases gephyrin phosphorylation at Ser270: a modification regulating postsynaptic GABAAR density. *Biol Chem.* doi: 10.1515/hsz-2021-0153.
5. Jeong, J. et al. (2021). Combination of cell signaling molecules can facilitate MYOD1-mediated myogenic transdifferentiation of pig fibroblasts. *J Anim Sci Biotechnol.* **12**(1):64. doi: 10.1186/s40104-021-00583-1.
6. Martin-Jaular, L. et al. (2021). Unbiased proteomic profiling of host cell extracellular vesicle composition and dynamics upon HIV-1 infection. *EMBO J.* doi: 10.15252/emboj.2020105492.
7. Xie, W. et al. (2021). Pterostilbene accelerates wound healing by modulating diabetes-induced estrogen receptor β suppression in hematopoietic stem cells. *Burns Trauma.* doi: 10.1093/burnst/tkaa045.
8. Sinha, A. et al. (2020). RAD51-Mediated DNA Homologous Recombination Is Independent of PTEN Mutational Status. *Cancers (Basel).* **12**(11):E3178. doi: 10.3390/cancers12113178.
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. Yoshihara, E. et al. (2020). Immune-evasive human islet-like organoids ameliorate diabetes. *Nature.* doi: 10.1038/s41586-020-2631-z.
11. Baci, D. et al. (2020). Extracellular Vesicles from Skeletal Muscle Cells Efficiently Promote Myogenesis in Induced Pluripotent Stem Cells. *Cells.* **9**(6):E1527. doi: 10.3390/cells9061527.
12. Smolková, K. et al. (2020). SIRT3 and GCN5L regulation of NADP⁺- and NADPH-driven reactions of mitochondrial isocitrate dehydrogenase IDH2. *Sci Rep.* **10**(1):8677. doi: 10.1038/s41598-020-65351-z.
13. Lucchini, F.C. et al. (2020). ASK1 inhibits browning of white adipose tissue in obesity. *Nat Commun.* **11**(1):1642. doi: 10.1038/s41467-020-15483-7.
14. Hsu, W.L. et al. (2020). Restoring Wnt6 signaling ameliorates behavioral deficits in MeCP2 T158A mouse model of Rett syndrome. *Sci Rep.* **10**(1):1074. doi: 10.1038/s41598-020-57745-w.
15. Harten, I.A. et al. (2020). The Synthesis and Secretion of Versican Isoform V3 by Mammalian Cells: A Role for N-linked Glycosylation. *Matrix Biol.* pii: S0945-053X(20)30008-1. doi: 10.1016/j.matbio.2020.01.002.
16. Yanai, M. et al. (2019). ADAR2 is involved in self and nonself recognition of Borna disease virus genomic RNA in the nucleus. *J Virol.* pii: JVI.01513-19. doi: 10.1128/JVI.01513-19.
17. Rubio, R.M. et al. (2019). Inhibition of ULK1 and Beclin1 by an α -herpesvirus Akt-like Ser/Thr kinase limits autophagy to stimulate virus replication. *Proc Natl Acad Sci U S A.* pii: 201915139. doi: 10.1073/pnas.1915139116.

18. Madan, E. et al. (2019). HIF-transcribed p53 chaperones HIF-1 α . *Nucleic Acids Res.* pii: gkz766. doi: 10.1093/nar/gkz766.
19. Zhou, Y. et al. (2019). Single-Cell Transcriptomic Analyses of Cell Fate Transitions during Human Cardiac Reprogramming. *Cell Stem Cell.* pii: S1934-5909(19)30224-3. doi: 10.1016/j.stem.2019.05.020.
20. Kobayashi, M. et al. (2019). AGO CLIP Reveals an Activated Network for Acute Regulation of Brain Glutamate Homeostasis in Ischemic Stroke. *Cell Rep.* **28**(4):979-991.e6. doi: 10.1016/j.celrep.2019.06.075.
21. Peng, D. et al. (2019). NRF2 antioxidant response protects against acidic bile salts-induced oxidative stress and DNA damage in esophageal cells. *Cancer Lett.* pii: S0304-3835(19)30326-X. doi: 10.1016/j.canlet.2019.05.031.
22. Lau, M.T. et al. (2019). Molecular dissection of box jellyfish venom cytotoxicity highlights an effective venom antidote. *Nat Commun.* **10**(1):1655. doi: 10.1038/s41467-019-09681-1.
23. Bhattacharya, S. et al. (2019). Lentiviral-mediated delivery of classical swine fever virus Erns gene into porcine kidney-15 cells for production of recombinant ELISA diagnostic antigen. *Mol Biol Rep.* doi: 10.1007/s11033-019-04829-0.
24. Navinés-Ferrer, A. et al. (2018). MRGPRX2-mediated mast cell response to drugs used in perioperative procedures and anaesthesia. *Sci Rep.* **8**(1):11628. doi: 10.1038/s41598-018-29965-8.
25. Meier, S. et al. (2018). Non-invasive detection of adeno-associated viral gene transfer using a genetically encoded CEST-MRI reporter gene in the murine heart. *Sci Rep.* **8**(1):4638. doi: 10.1038/s41598-018-22993-4.
26. Ghosh, N. & Banerjee, E.R. (2017). A Review on Various Tissue Engineering Techniques to Induce Differentiation of Pluripotent Stem Cells. *Medical Glory.* **1**(2):130-149.
27. Agudelo, D. et al. (2017). Marker-free coselection for CRISPR-driven genome editing in human cells. *Nat Methods.* **14**(6):615–620. doi:10.1038/nmeth.4265.
28. Zheng, L. et al. (2017). Lym-1 Chimeric Antigen Receptor T Cells Exhibit Potent Anti-Tumor Effects against B-Cell Lymphoma. *Int J Mol Sci.* **18**(12). pii: E2773. doi: 10.3390/ijms18122773.
29. Hsu, W.L. et al. (2017). Smad4 SUMOylation is essential for memory formation through upregulation of the skeletal myopathy gene TPM2. *BMC Biol.* **15**(1):112. doi: 10.1186/s12915-017-0452-9.
30. Tao, C.C. et al. (2017). Epigenetic regulation of HDAC1 SUMOylation as an endogenous neuroprotection against A β toxicity in a mouse model of Alzheimer's disease. *Cell Death Differ.* doi: 10.1038/cdd.2016.161.

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

©2006-2021: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.