

# Epi5™ Episomal iPSC Reprogramming Kit

## USER GUIDE

For efficient, integration-free reprogramming of somatic cells into induced pluripotent stem cells (iPSC)

Catalog Number A15960

Publication Number MAN0008352

Revision A.0

Information in this guide is subject to change without notice.

## DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

## Revision History

Revision	Date	Description
A.0	March 2016	Rebrand; update legal boilerplates; update Methods
1.0	11 June 2013	Baseline for this revision

## LIMITED USE LABEL LICENSE No. 465: Epi5™ Vectors

**Notice to Purchaser:** The transfer of this product conveys to the recipient user the limited, non-transferable right to use the transferred amount of product to perform internal research for the sole benefit of the recipient user and for educational purposes. No right to resell this product or any of its components, or iPS cells generated by use of the product, or modifications (progenitors or fully differentiated cells) thereof is conveyed expressly, by implication, or by estoppels. This product, iPS cells made using this product and their modifications shall not be administered to – (a) human subjects, including for human clinical use and/or (b) animals for veterinary use (i.e., not for research) – for therapeutic, diagnostic or prophylactic purposes including but not limited to clinical applications, cell therapy, transplantation and/or regenerative medicine, nor shall be used for the creation of human embryos, and/or admixed embryos with embryos of animals including humans for any purpose including for research. This product, iPS cells made using this product and their modifications is for internal research purposes and is not for use in any of the following commercial applications: (i) use of iPS cells and progeny (but not modifications) for manufacture or quality control of any product; (ii) use of cells to provide a service, information or data to a third party for financial gain (financial benefit exceeds the cost of operating the transaction), except when performed by a not-for-profit research entity (not affiliated with a for-profit organization) not for financial gain, or performed on behalf of a third party having an appropriate license\*; (iii) use of cells by or on behalf of recipient user for screening or later stage development of therapeutics, diagnostics, prophylactics (e.g., hit-to-lead, lead optimization), except when performed by or on behalf of a not-for-profit research entity for internal research and not for financial gain, or when performed in accordance with a separate Limited Sub License Agreement for use of LTC Pluripotent Cells executed in writing between recipient user and LTC; (iv) sale, lease distribution, transfer or license of cells to third parties for financial gain; (v) any sale, lease, distribution, transfer or license of iPS cells and progeny (but not modifications) by recipient user to for-profit entities, even if for research use. Any person or entity that wishes to perform commercial applications described above using this product, iPS cells derived from the use of this product, or modifications thereof, is required to take a license to cover the appropriate commercial application. An entity that wants to transfer iPS cells or progeny to a for-profit recipient shall notify the for-profit recipient that such recipient is required to take the appropriate license to use the materials for internal research of any type.

Notwithstanding the foregoing, the following activities are not considered commercial applications for purposes of this label license:

- (a) basic research, including, without limitation, target discovery, target validation and assay development;
- (b) transfer of cells to not-for-profit research entity for its internal research not for financial gain;
- (c) compound screening and safety testing for development of therapeutics, diagnostics and prophylactics by academic and not-for-profit research entities for their non-commercial internal research;
- (d) license or commercialization of research results except where such results are drugs or drug candidates, iPS cells or modifications, or where such license or commercialization uses iPS cells or progeny.

Other than rights granted herein, no other right, express or implied, is conveyed by the sale of this product.

\*For information on obtaining additional rights, please contact [outlicensing@lifetechnologies.com](mailto:outlicensing@lifetechnologies.com) or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, CA 92008.

Notice: If iPS cells and derivatives (hereafter “Materials”) generated by use of this product are transferred to other users in accordance with the terms of the label license for reprogramming technology set forth above and accompanying the product (hereafter “Label License”), the transferring party should notify the recipients of such Materials of these terms by transferring a copy of the Label License to the recipients.

## Corporate entity

Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

## Trademarks

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Parafilm is a registered trademark of Bemis Company, Inc. B-27 is a registered trademark of the Board of Trustees of Southern Illinois University.

© 2016 Thermo Fisher Scientific. All rights reserved.

# Contents

<b>Product information</b> .....	<b>4</b>
Product description.....	4
Contents and storage .....	4
Description of the system .....	5
Materials and equipment required .....	7
<b>Methods</b> .....	<b>9</b>
Guidelines for generating iPSCs.....	9
Reprogram fibroblasts .....	10
Reprogram StemPro™ CD34 <sup>+</sup> Cells (Feeder-Free Culture).....	13
Reprogramming StemPro™ CD34 <sup>+</sup> Cells (Feeder-Dependent Culture) .....	16
Identify iPSC colonies .....	19
Live stain with antibodies .....	21
Pick iPSC colonies .....	22
Detect episomal vectors by PCR.....	24
<b>Appendix A: Recipes</b> .....	<b>27</b>
Media and reagents.....	27
<b>Appendix B: Prepare culture vessels</b> .....	<b>30</b>
Coat culture vessels with Geltrex™ matrix.....	30
MEF culture dishes.....	31
<b>Appendix C: Cryopreserve iPSCs</b> .....	<b>33</b>
Freeze iPSCs (Feeder-Free Culture) .....	33
Freeze iPSCs (Feeder-Dependent Culture) .....	34
<b>Appendix D: Ordering information</b> .....	<b>36</b>
Accessory products .....	36
<b>Documentation and support</b> .....	<b>39</b>
Customer and technical support .....	39

# Product information

## Product description

---

### Induced pluripotent stem cells (iPSC)

Induced pluripotent stem cells (iPSCs) are genetically reprogrammed adult cells which exhibit a pluripotent stem cell-like state similar to embryonic stem cells.

The Invitrogen™ Epi5™ Episomal iPSC Reprogramming Kit contains an optimized mixture of vectors that can reprogram somatic cells to iPSCs without integration. The Epi5™ Episomal iPSC Reprogramming Kit contains **five** reprogramming vectors in two separate tubes. Tube A contains an optimized mixture of three episomal vectors that deliver the five reprogramming factors Oct4, Sox2, Klf4, L-Myc, and Lin28. Tube B contains an optimized mixture of two vectors expressing mp53DD (a dominant negative mutation of p53) and EBNA1, which together improve the reprogramming efficiency of the system.

The Epi5™ Episomal iPSC Reprogramming Kit contains sufficient material for 20 reprogramming experiments.

## Contents and storage

---

### Epi5™ Episomal iPSC Reprogramming Kit (Cat. No. A15960)

Contents	Amount	Vectors	Genes	Storage
Epi5™ Reprogramming Vectors (Tube A)	20 µL	pCE-hOCT3/4	Oct4	-20°C
		pCE-hSK	Sox2, Klf4	
		pCE-hUL	L-Myc, Lin28	
Epi5™ p53 & EBNA Vectors (Tube B)	20 µL	pCE-mP53DD	mp53DD	
		pCXB-EBNA1	EBNA1	

## Description of the system

---

<b>Induced pluripotent stem cells (iPSC)</b>	<p>Induced pluripotent stem cells (iPSCs) are genetically reprogrammed adult cells which exhibit a pluripotent stem cell-like state similar to embryonic stem cells. While these artificially generated cells are not known to exist in the human body, they show qualities remarkably similar to those of embryonic stem cells (ESC); thus, they are an invaluable new source of pluripotent cells for drug discovery, cell therapy, and basic research.</p> <p>There are multiple methods to generate iPSCs, including retrovirus-mediated gene transfection and chemical induction. While retroviral vectors require integration into host chromosomes to express reprogramming genes, DNA-based vectors and plasmid vectors exist episomally and do not require integration. The Epi5™ Episomal iPSC Reprogramming Kit contains an optimized mixture of <b>five</b> vectors that can reprogram somatic cells to iPSCs without integration.</p>
<b>Epi5™ Reprogramming Vectors</b>	<p>The Epi5™ Reprogramming Vectors contain an optimized mixture of three episomal vectors with an oriP/EBNA-1 (Epstein-Barr nuclear antigen-1) backbone for delivering the reprogramming genes, Oct4, Sox2, Lin28, L-Myc, and Klf4. High transfection efficiency due to oriP/EBNA-1 mediated nuclear import and retention of vector DNA allows iPSC derivation in a single transfection. In addition, silencing of the viral promoter driving EBNA-1 expression and the loss of the episomes at a rate of ~5% per cell cycle due to defects in vector synthesis and partitioning allows the removal of episomal vectors from the iPSCs without any additional manipulation.</p>
<b>Epi5™ p53 &amp; EBNA Vectors</b>	<p>Epi5™ p53 &amp; EBNA Vectors provide additional improvements to the reprogramming system.</p> <p>The p53 protein is highly involved in cell cycle regulation and tumor suppression. With the activation of cytotoxic responses, p53 expression results in cell cycle arrest or cell death to prevent further complications within the system. Knockdown of p53 has been shown to improve reprogramming efficiencies as well as to prevent differentiation via the introduction of a variety of knockdown agents. The mp53DD is a dominant negative mutation of the p53 protein providing higher efficiency knockdown than traditional shRNA systems. The presence of this gene in an episomal system allows for transient expression of the dominant negative mutant over an extended period of time, diluting out with normal passaging of iPSC.</p> <p>Supplemental expression of the EBNA1 gene from a vector solely dedicated to this purpose allows for high expression of plasmids containing the origin of replication present on the reprogramming plasmids. This plasmid, however, lacks the OriP gene, thus providing only transient expression for the early stages of reprogramming.</p>

## Comparison of episomal reprogramming vectors

The table below provides a comparison of the Episomal iPSC Reprogramming Vectors (Cat. no. A14703) and the Epi5™ Episomal iPSC Reprogramming Kit (Cat. no. A15960).

	Episomal iPSC Reprogramming Vectors	Epi5™ Episomal iPSC Reprogramming Kit
<b>Genes</b>	Oct4, Sox2, Nanog, Lin28, Klf4, and L-Myc	Oct4, Sox2, Lin28, L-Myc, Klf4, mp53DD, additional EBNA
<b>Vector size</b>	12–17.5 kb	6–11 kb
<b>Polycistronic elements</b>	IRES element	2A sequence
<b>Promoter</b>	EF1 $\alpha$	CAG
<b>Small molecules</b>	required	not required
<b>Number of vectors</b>	3	5

## Advantages of Epi5™ Episomal iPSC Reprogramming Kit

- As a transgene-free and viral-free reprogramming system, it is a safe alternative to other reprogramming methods such as lentiviral delivery for all stages of iPSC research.
- Allows the reprogramming of a variety of somatic cell types and provides flexibility in somatic cell selection.
- Optimized for feeder-free reprogramming, enabling defined and feeder-free reprogramming when used with the Essential 8™ Medium.
- High transfection efficiency due to oriP/EBNA-1 mediated nuclear import and retention of vector DNA allows iPSC derivation in a single transfection.
- The episomal vectors are removed from the iPSCs without any further manipulation due to the silencing of the viral promoter driving EBNA-1 expression and the loss of the episomes due to defects in vector synthesis and partitioning.
- Do not require the use of small molecules for reprogramming.

## Materials and equipment required

---

- Cells and vectors**
- Epi5™ Episomal iPSC Reprogramming Kit (Cat. No. A15960)
  - Human neonatal foreskin fibroblast cells (strain BJ; ATCC no. CRL2522)
- Media and reagents**
- Dulbecco's Modified Eagle Medium (DMEM) with GlutaMAX™-I (high glucose) (Cat. No. 10569)
  - KnockOut™ DMEM/F-12 (Cat. No. 12660-012)
  - Fetal Bovine Serum (FBS), ESC-Qualified (Cat. No. 16141-079)
  - MEM Non-essential Amino Acids (NEAA) (Cat. No. 11140-050)
  - Basic Fibroblast Growth Factor (bFGF), recombinant human (Cat. No. PHG0264)
  - Bovine Albumin Fraction V Solution (BSA) (Cat. No. 15260-037)
  - Essential 8™ Medium, consisting of Essential 8™ Basal Medium and Essential 8™ Supplement (50X) (Cat. No. A1517001)
  - DMEM/F-12 with HEPES (Cat. No. 11330-057)
  - N-2 Supplement (100X) (Cat. No. 17502-048)
  - B-27™ Serum-Free Supplement (50X) (Cat. No. 17504-044)
  - GlutaMAX™-I Supplement (Cat. No. 35050-061)
  - β-mercaptoethanol (Cat. No. 21985-023)
  - Geltrex™ LDEV-Free hESC-Qualified Reduced Growth Factor Basement Membrane Matrix (Cat. No. A1413301)
  - 0.05% Trypsin-EDTA (1X), Phenol Red (Cat. No. 25300-054)
  - UltraPure™ 0.5 M EDTA, pH 8.0 (Cat. No. 15575-020)
  - Dulbecco's PBS (DPBS) without Calcium and Magnesium (Cat. No. 14190144)
- Characterization reagents**
- Use one of the following:
- TRA-1-60 Alexa Fluor™ 488 Conjugate Kit for Live Cell Imaging (Cat. No. A25618)
  - TRA-1-60 Alexa Fluor™ 594 Conjugate Kit for Live Cell Imaging (Cat. No. A24882)
  - TRA-1-60 Alexa Fluor™ 555 Conjugate Kit for Live Cell Imaging (Cat. No. A24879)
- Reagents for PCR detection of episomal vectors**
- CellsDirect Resuspension & Lysis Buffers (Cat. No. 11739-010)
  - AccuPrime™ Taq DNA Polymerase High Fidelity (Cat. No. 12346-086)
  - Forward and Reverse primers for PCR (see page 26)

## Equipment

- Electroporation instrument.  
**Note:** The protocols in this user guide describe electroporation using the Neon™ Transfection System (Cat. no. MPK5000) and the Neon™ Transfection System 10 µL Tip Kit (Cat. no. MPK1025). Electroporation parameters must be optimized for other electroporation devices.
- Sterile cell culture hood (i.e., biosafety cabinet) equipped with a stereomicroscope
- Inverted microscope
- Incubator set at 37°C, 5% CO<sub>2</sub>
- Water bath set at 37°C
- Sterile serological pipettes (5-mL, 10-mL)
- Centrifuge
- 15-mL centrifuge tubes
- 60-mm and 100-mm tissue culture-treated dishes
- 6-well tissue culture-treated plates
- 25 gauge 1 inch needle



## Guidelines for generating iPSCs

---

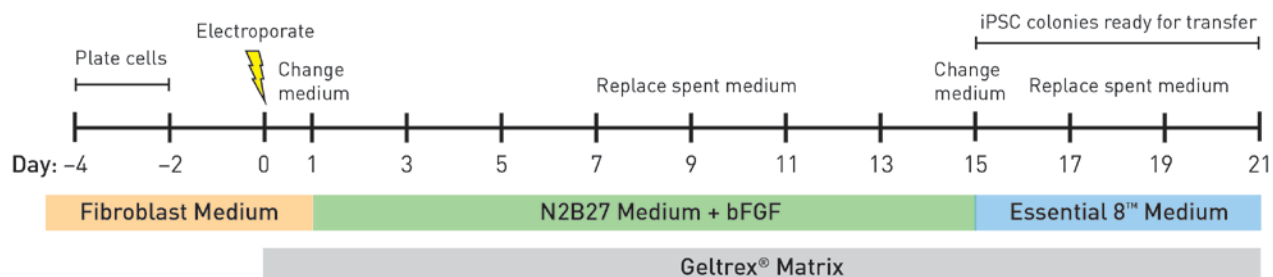
### Experiment guidelines

- To maintain sterile culture conditions, carry out all of the procedures in this manual using sterile laboratory practices in a laminar flow hood.
- You can use the Epi5™ Episomal iPSC Reprogramming Kit to reprogram a wide range of cell types in proliferative and quiescent states. However, the reprogramming efficiency may vary between different cell types.
- Each reprogramming experiment requires  $1 \times 10^5$  cells and 1  $\mu$ L each of the Epi5™ Reprogramming Vectors and the Epi5™ p53 & EBNA Vectors.
- The Epi5™ Episomal iPSC Reprogramming Kit contains sufficient material for 20 reprogramming experiments. Before use, briefly centrifuge the thawed vectors to collect them at the bottom of the tube.
- Cells should be approximately 75–90% confluent on the day of transfection (Day 0). Since overconfluency results in decreased transfection efficiency, we recommend replating the cells to achieve 75–90% confluency, if they have become overconfluent during culturing.
- We recommend reprogramming patient cells at the earliest passage possible. However, it is important to have the cells growing and healthy, which can take between 1–4 weeks. The cells are usually ready to reprogram once they have gone through a total of 3–4 passages.
- Prior to starting your experiments, ensure that the media are equilibrated to 37°C and appropriately gassed.
- You do not need to manipulate the culture conditions to remove the reprogramming vectors. In the absence of drug selection, approximately 5% of the episomal vectors are lost each cell cycle due to silencing of the viral promoter driving EBNA-1 expression and defects in vector synthesis and partitioning.
- Newly derived iPSC lines may contain a fair amount of differentiation through passage 4. It is not necessary to remove differentiated material prior to passaging. By propagating/splitting the cells the overall culture health should improve throughout the early passages.

# Reprogram fibroblasts

## Workflow

The major steps required for reprogramming human neonatal foreskin fibroblast cells to generate iPSCs using the Epi5™ Episomal iPSC Reprogramming Kit are shown below. Note that the timeline is provided as a guideline for experimental planning; actual timeline can vary based on the cell type and experimental conditions.



## Fibroblast reprogramming timeline

**Day -4 to -2:** Plate human fibroblasts into a T75 flask in Fibroblast Medium so that they are 75–90% confluent on the day of transfection (Day 0).

**Day 0:** Transfect the cells using the Neon™ Transfection System. Plate transfected cells onto Geltrex™ matrix-coated culture dishes, and incubate them overnight in complete Fibroblast Medium.

**Day 1 to 14:** Change the medium to N2B27 Medium supplemented with 100 ng/mL bFGF; replace the spent medium every other day.

**Day 15:** Change the medium to Essential 8™ Medium and monitor the culture vessels for the emergence of iPSC colonies.

**Day 21:** Pick and transfer undifferentiated iPSCs onto fresh Geltrex™ matrix-coated culture dishes for expansion.

## Media for reprogramming fibroblasts

For optimal reprogramming efficiency with the Epi5™ Episomal iPSC Reprogramming Kit, culture the cells in Fibroblast Medium on Geltrex™ matrix-coated culture vessels until the day of transfection.

After transfection, allow the cells to recover in Fibroblast Medium overnight, and then switch to N2B27 Medium supplemented with 100 ng/mL bFGF.

After 15 days of culture in Fibroblast Medium and on Geltrex™ matrix, transition the reprogrammed cells into Essential 8™ Medium, a serum-free, xeno-free medium, and continue culturing them on Geltrex™ matrix-coated culture vessels until iPSCs emerge and are ready to be picked.

See page 27 for instructions on preparing media and reagents.

## Fibroblast reprogramming protocol

The following reprogramming protocol has been optimized for human neonatal foreskin fibroblast cells (strain BJ; ATCC no. CRL2522). We recommend that you optimize the protocol for your cell type.

### Day -4 to -2: Seed cells

2–4 days before transfection, plate human fibroblast cells in Fibroblast Medium (see page 27 for recipe) into a T75 flask. Cells should be approximately 75–90% confluent on the day of transfection (Day 0).

**Note:** Growth rate is dependent on the cell line and culture conditions. Depending on the seeding density and culture conditions, the cells may take up to 5 days to reach 75–90% confluency.

**Note:** Because overconfluency results in decreased transfection efficiency, we recommend

replating your cells to achieve 75–90% confluency if your cells have become overconfluent during culturing.

**Day 0:  
Prepare the cells  
for transfection**

Gentle handling of the cells prior to transfection is essential for the success of the transfection procedure.

- Aspirate the medium from Geltrex™ matrix-coated 6-well plates (see page 30) and replace with 2 mL of fresh Fibroblast Medium per well. Place the coated plates at 37°C until ready for use.  
**Note:** One Geltrex™ matrix-coated 6-well plate can be used for six transfections.
- Thaw the Epi5™ vectors at 37°C and place them on wet ice until ready for use. Before use, briefly centrifuge the thawed vectors to collect them at the bottom of the tube.
- Wash the cells in DPBS without calcium and magnesium.
- Add 3 mL of 0.05% Trypsin/EDTA to the culture flask containing the cells.
- Incubate the flask at 37°C for approximately 4 minutes.
- Add 6 mL of Fibroblast Medium to each flask. Tap the plate against your hand to ensure cells have been dislodged from the flask, and carefully transfer cells into an empty, sterile 15-mL conical tube.
- Wash the culture flask with an additional 4 mL of Fibroblast Medium and add to the wash to the same 15-mL conical tube to ensure that all of the cells have been transferred.
- Remove a 20-µL sample to perform a viable cell count and calculate the number of transfections to be performed. You will need  $1 \times 10^5$  cells per transfection.
- Centrifuge the cells at  $200 \times g$  for 5 minutes at room temperature.
- Carefully aspirate most of the supernatant, leaving approximately 100–200 µL behind. Remove the remaining medium with a 200-µL pipette.

**Day 0:  
Transfect cells  
using the Neon™  
Transfection System**

- Resuspend the cell pellet in Resuspension Buffer R (included with Neon™ Transfection kits) at a final concentration of  $1.0 \times 10^7$ – $1.4 \times 10^7$  cells/mL.
- Transfer the cells (10 µL per transfection reaction) to a sterile 1.5-mL microcentrifuge tube.
- Turn on the Neon™ unit and enter the following electroporation parameters in the Input window.

Pulse voltage	Pulse width	Pulse number	Cell density	Tip type
1650 V	10 ms	3	$1 \times 10^7$ cells/mL	10-µL

- Fill the Neon™ Tube with 3 mL Electrolytic Buffer (use Buffer E for the 10-µL Neon™ Tip).
- Insert the Neon™ Tube into the Neon™ Pipette Station until you hear a click.
- Transfer 1 µL of Epi5™ Reprogramming Vectors to the tube containing cells.
- Transfer 1 µL of Epi5™ p53 & EBNA Vectors to the tube containing cells and mix gently.
- Insert a 10-µL Neon™ Tip into the Neon™ Pipette.
- Press the push-button on the Neon™ Pipette to the first stop and immerse the Neon™ Tip into the cell-DNA mixture. Slowly release the push-button on the pipette to aspirate the cell-DNA mixture into the Neon™ Tip.

**Note:** Avoid air bubbles during pipetting to avoid arcing during electroporation. If you notice air bubbles in the tip, discard the sample and carefully aspirate fresh

sample into the tip again without any air bubbles.

10. Insert the Neon™ Pipette with the sample vertically into the Neon™ Tube placed in the Neon™ Pipette Station until you hear a click.
11. Ensure that you have entered the appropriate electroporation parameters and press Start on the Neon™ touchscreen to deliver the electric pulse.  
**Note:** After the electric pulse is delivered, the touchscreen displays “Complete” to indicate that electroporation is complete.
12. Remove the Neon™ Pipette from the Neon™ Pipette Station and immediately transfer the samples from the Neon™ Tip into one well of the pre-warmed Geltrex™ matrix-coated 6-well plate containing Fibroblast Medium.  
**Note:** Evenly distribute cells over the well in a dropwise manner.
13. Discard the Neon™ Tip into an appropriate biological hazardous waste container.
14. Repeat the process for any additional samples. Do not use Neon™ tip more than twice.
15. Incubate the plates at 37°C in a humidified CO<sub>2</sub> incubator overnight.

**Day 1:**

**Switch to N2B27 Medium**

1. Aspirate the spent Fibroblast Medium from the plates.
2. Add 2 mL of N2B27 Medium supplemented with bFGF (added freshly prior to use) to each well (see page 28 for recipe).
3. Replace the spent medium every other day, up to day 15 post-transfection.

**Day 15:**

**Switch to**

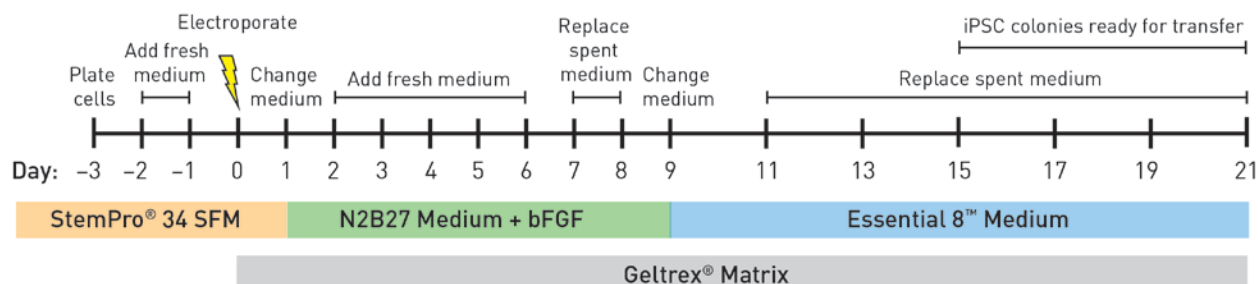
**Essential 8™ Medium**

1. Aspirate the spent medium and replace with Essential 8™ Medium (see page 28 for recipe). Resume medium changes every day.
2. Observe the plates every other day under a microscope for the emergence of cell clumps indicative of transformed cells (see **Identify iPSC colonies**, page 19). Within 15 to 21 days of transfection, the iPSC colonies will grow to an appropriate size for transfer.

# Reprogram StemPro™ CD34<sup>+</sup> Cells (Feeder-Free Culture)

## Workflow

The major steps required for reprogramming StemPro™ CD34<sup>+</sup> cells to generate iPSCs using the Epi5™ Episomal iPSC Reprogramming Kit and the subsequent culture of the reprogrammed cells under feeder-free conditions are shown below. Note that the timeline is provided as a guideline for experimental planning; actual timeline can vary based on the cell type and experimental conditions.



## StemPro™ CD34<sup>+</sup> (feeder-free) reprogramming timeline

**Day -3:** Plate StemPro™ CD34<sup>+</sup> cells into a 24-well plate in complete StemPro™-34 medium containing cytokines (i.e., SCF, IL-3, and GM-CSF).

**Day 0:** Transfect the cells using the Neon™ Transfection System. Plate transfected cells onto Geltrex™ matrix-coated culture dishes, and incubate them overnight in complete StemPro™-34 medium containing cytokines.

**Day 1:** Change the medium to N2B27 Medium supplemented with 100 ng/mL bFGF; replace the spent medium every other day.

**Day 2 to 6:** Replenish the culture every day by adding fresh medium.

**Day 7 to 8:** Replace the spent medium with fresh medium every day.

**Day 9:** Change the medium to Essential 8™ Medium and monitor the culture vessels for the emergence of iPSC colonies. Replace the spent medium every other day.

**Day 15 to 21:** Pick and transfer undifferentiated iPSCs onto fresh Geltrex™ matrix-coated culture dishes for expansion.

## Media for reprogramming StemPro™ CD34<sup>+</sup> cells (feeder-free)

For optimal reprogramming efficiency with the Epi5™ Episomal iPSC Reprogramming Kit, culture the cells in complete StemPro™-34 medium containing cytokines (i.e., SCF, IL-3, and GM-CSF) until the day of transfection. After transfection, allow the cells to recover in complete StemPro™-34 medium with cytokines on Geltrex™ matrix-coated culture vessels overnight, and then switch to N2B27 Medium supplemented with 100 ng/mL bFGF.

After 9 days of culture in supplemented N2B27 Medium on Geltrex™ matrix-coated culture vessels, transition the reprogrammed cells into Essential 8™ Medium, and continue culturing them on Geltrex™ matrix-coated culture vessels until iPSCs emerge and are ready to be picked.

See page 27 for instructions on preparing media and reagents.

## StemPro™ CD34<sup>+</sup> reprogramming protocol (feeder-free culture)

The following reprogramming protocol has been optimized for StemPro™ CD34<sup>+</sup> cells (Cat. no. A14059), which are human hemopoietic progenitor cells (HPCs) derived from the human umbilical cord blood of mixed donors. We recommend that you optimize the protocol for your cell type.

**Day –3 to –1:  
Seed and expand  
cells**

- 3 days before transduction (day –3), thaw 1 vial of StemPro™ CD34<sup>+</sup> cells (0.5 × 10<sup>6</sup> cells) and gently transfer into a 15-mL sterile, conical tube. Drop-wise add 10 mL of StemPro™-34 SFM to the cells while gently agitating to mix.
- Centrifuge the cell suspension at 200 × g for 10 minutes, discard the supernatant, and resuspend the cells in 1 mL of complete StemPro™-34 medium containing cytokines (i.e., SCF, IL-3, and GM-CSF) into one well of a 24-well plate (see page 28).  
**Note:** The recommended final concentrations for the cytokines are 100 ng/mL for SCF, 50 ng/mL for IL-3, and 25 ng/mL for GM-CSF.  
**Note:** We recommend using the wells in the middle section of the 24-well plate to prevent excessive evaporation of the medium during incubation.
- 2 days before transduction (day –2), add 0.5 mL of fresh complete StemPro™-34 medium containing cytokines without disturbing the cells.
- One day before transduction (day –1), gently remove 0.5 mL of the medium, and add 1.0 mL of fresh complete StemPro™-34 medium containing cytokines without disturbing the cells.

**Day 0:  
Transfect cells  
using the Neon™  
Transfection  
System**

Gentle handling of the cells prior to transfection is essential for the success of the transfection procedure.

- Aspirate the medium from Geltrex™ matrix-coated 6-well plates (see page 30) and replace with 2 mL of complete StemPro™-34 medium containing cytokines (see page 28) per well. Place the coated plates at 37°C until ready for use.  
**Note:** One Geltrex™ matrix-coated 6-well plate can be used for six transfections.
- Thaw the Epi5™ vectors at 37°C and place them on wet ice until ready for use. Before use, briefly centrifuge the thawed vectors to collect them at the bottom of the tube.
- Count the number of cells using a hemacytometer or the Countess™ Automated Cell Counter to determine the viability and total number of cells; the cells should have more than doubled in number.
- Carefully aspirate most of the supernatant, leaving approximately 100–200 µL behind. Remove the remaining medium with a 200-µL pipette.
- Resuspend the cell pellet in Resuspension Buffer T (included with Neon™ Transfection kits) at a final concentration of 1.0 × 10<sup>7</sup> –1.4 × 10<sup>7</sup> cells/mL.  
**Note:** Buffer T is required for blood cell transfections.

**Day 0:  
Transfect cells  
using the Neon™  
Transfection  
System, continued**

- Transfer the cells (10 µL per transfection reaction) to a sterile 1.5-mL microcentrifuge tube.
- Turn on the Neon™ unit and enter the following electroporation parameters in the Input window.

Pulse voltage	Pulse width	Pulse number	Cell density	Tip type
1650 V	10 ms	3	1 × 10 <sup>7</sup> cells/mL	10-µL

- Fill the Neon™ Tube with 3 mL Electrolytic Buffer (use Buffer E for the 10-µL Neon™ Tip).
- Insert the Neon™ Tube into the Neon™ Pipette Station until you hear a click.
- Transfer 1 µL of Epi5™ Reprogramming Vectors to the tube containing cells.
- Transfer 1 µL of Epi5™ p53 & EBNA Vectors to the tube containing cells and mix gently.

12. Insert a 10- $\mu$ L Neon™ Tip into the Neon™ Pipette.
13. Press the push-button on the Neon™ Pipette to the first stop and immerse the Neon™ Tip into the cell-DNA mixture. Slowly release the push-button on the pipette to aspirate the cell-DNA mixture into the Neon™ Tip.
 

**Note:** Avoid air bubbles during pipetting to avoid arcing during electroporation. If you notice air bubbles in the tip, discard the sample and carefully aspirate fresh sample into the tip again without any air bubbles.
14. Insert the Neon™ Pipette with the sample vertically into the Neon™ Tube placed in the Neon™ Pipette Station until you hear a click.
15. Ensure that you have entered the appropriate electroporation parameters and press **Start** on the Neon™ touchscreen to deliver the electric pulse.
 

**Note:** After the electric pulse is delivered, the touchscreen displays “Complete” to indicate that electroporation is complete.
16. Remove the Neon™ Pipette from the Neon™ Pipette Station and immediately transfer the samples from the Neon™ Tip into one well of the pre-warmed Geltrex™ matrix-coated 6-well plate containing 2 mL of complete StemPro™-34 medium with cytokines.
 

**Note:** Evenly distribute cells over the well in a dropwise manner.
17. Discard the Neon™ Tip into an appropriate biological hazardous waste container.
18. Repeat the process for any additional samples. Do not use Neon™ tip more than twice.
19. Incubate the plates at 37°C in a humidified CO<sub>2</sub> incubator overnight.

**Day 1:  
Switch to N2B27  
Medium**

1. Do not aspirate the spent medium from the plates. Add 1 mL of N2B27 Medium supplemented with 100 ng/mL bFGF (added freshly prior to use; see page 28 for recipe) to each well.

**Day 2 to 6:  
Add fresh N2B27  
Medium**

1. Carefully aspirate 0.5–1 mL of the supernatant from the top of the well, trying to not remove any of the suspension StemPro™ CD34<sup>+</sup> cells while they are still attaching.
2. Add 0.5–1 mL of N2B27 medium supplemented with 100 ng/mL bFGF every day, up to day 6 post-transfection.

**Day 7 to 9:  
Replace spent  
N2B27 Medium**

1. Carefully aspirate the spent medium and replace with 2 mL of N2B27 medium supplemented with 100 ng/mL bFGF.
2. Replace spent medium with 2 mL fresh medium every day, up to day 9 post-transfection.

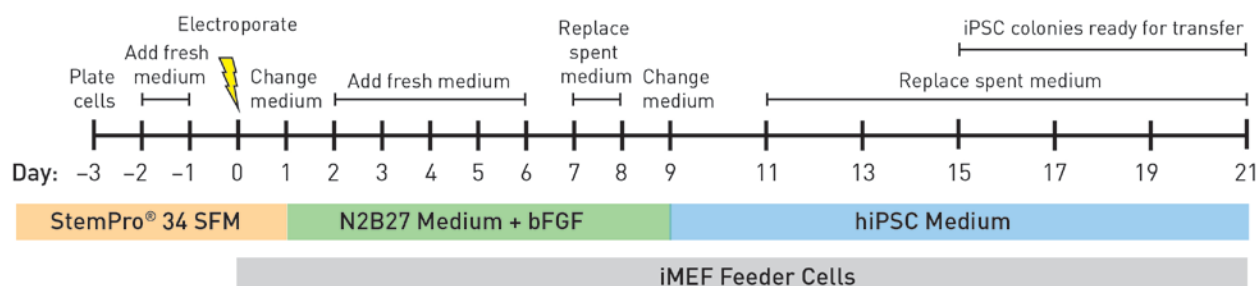
**Day 9:  
Switch to  
Essential 8™  
Medium**

1. Aspirate the spent N2B27 medium and replace it with complete Essential 8™ Medium (see page 28 for recipe). Resume medium changes every day.
2. Observe the plates every other day under a microscope for the emergence of cell clumps indicative of transformed cells (see **Identify iPSC colonies**, page 19). Within 15 to 21 days of transfection, the iPSC colonies will grow to an appropriate size for transfer.

# Reprogramming StemPro™ CD34+ Cells (Feeder-Dependent Culture)

## Workflow

The major steps required for reprogramming StemPro™ CD34+ cells to generate iPSCs using the Epi5™ Episomal iPSC Reprogramming Kit and the subsequent culture of the reprogrammed cells on MEF feeder layers are shown below. Note that the timeline is provided as a guideline for experimental planning; actual timeline can vary based on the cell type and experimental conditions.



## StemPro™ CD34+ (feeder-dependent) reprogramming timeline

**Day -3:** Plate the StemPro™ CD34+ cells into a 24-well plate in complete StemPro™-34 medium containing cytokines (i.e., SCF, IL-3, and GM-CSF).

**Day 0:** Transfect the cells using the Neon™ Transfection System. Plate transfected cells onto MEF feeders, and incubate them overnight in complete StemPro™-34 medium containing cytokines.

**Day 1:** Change the medium to N2B27 Medium supplemented with 100 ng/mL bFGF.

**Day 2 to 6:** Replenish the culture every day by adding fresh medium.

**Day 7 to 8:** Replace the spent medium with fresh medium every day.

**Day 9:** Change the medium to human iPSC Medium and continue culturing the reprogrammed cells on MEF feeders, monitoring the culture for the emergence of iPSC colonies and replacing the spent medium every day.

**Day 15 to 21:** Pick and transfer undifferentiated iPSCs onto fresh culture dishes with MEF feeder layers for expansion.

## Media for reprogramming StemPro™ CD34+ cells (feeder-dependent)

For optimal reprogramming efficiency using the Epi5™ Episomal iPSC Reprogramming Kit, culture the cells in complete StemPro™-34 medium containing cytokines (i.e., SCF, IL-3, and GM-CSF) until the day of transfection.

After transfection, allow the cells to recover in complete StemPro™-34 medium with cytokines on MEF feeders overnight, and then switch to N2B27 Medium supplemented with 100 ng/mL bFGF.

After 9 days of culture in supplemented N2B27 Medium on MEF feeder layers, transition the reprogrammed cells into human iPSC Medium, and continue culturing them on MEF feeders until iPSCs emerge and are ready to be picked.

See page 27 for instructions on preparing media and reagents.

## StemPro™ CD34+ Reprogramming Protocol (Feeder-Dependent Culture)

The following reprogramming protocol has been optimized for StemPro™ CD34+ cells (Cat. no. A14059), which are human hemopoietic progenitor cells (HPCs) derived from the human umbilical cord blood of mixed donors. We recommend that you optimize the protocol for your cell type.



**Day –3 to –1:  
Seed and expand  
cells**

- 3 days before transduction (day –3), thaw 1 vial of StemPro™ CD34+ cells ( $0.5 \times 10^6$  cells) and gently transfer into a 15-mL sterile, conical tube. Drop-wise add 1 mL of StemPro™-34 SFM to the cells while gently agitating to mix.
- Centrifuge the cell suspension at  $200 \times g$  for 10 minutes, discard the supernatant, and resuspend the cells in 1 mL of complete StemPro™-34 medium containing cytokines (i.e., SCF, IL-3, and GM-CSF) into one well of a 24-well plate (see page 28).  
**Note:** The recommended final concentrations for the cytokines are 100 ng/mL for SCF, 50 ng/mL for IL-3, and 25 ng/mL for GM-CSF.  
**Note:** We recommend using the wells in the middle section of the 24-well plate to prevent excessive evaporation of the medium during incubation.
- 2 days before transduction (day –2), add 0.5 mL of fresh complete StemPro™-34 medium containing cytokines without disturbing the cells.
- One day before transduction (day –1), gently remove 0.5 mL of the medium, and add 1.0 mL of fresh complete StemPro™-34 medium containing cytokines without disturbing the cells.

**Day 0:  
Transfect cells  
using the Neon™  
Transfection  
System**

Gentle handling of the cells prior to transfection is essential for the success of the transfection procedure.

- Aspirate the medium from the 6-well plates containing MEF feeder cells (see page 31) and replace it with 2 mL per well of complete StemPro™-34 medium containing cytokines (see page 28). Place the coated plates at 37°C until ready for use.  
**Note:** One 6-well plate with MEF feeder layers can be used for six transfections.
- Thaw the Epi5™ vectors at 37°C and place them on wet ice until ready for use. Before use, briefly centrifuge the thawed vectors to collect them at the bottom of the tube.
- Count the number of cells using a hemacytometer or the Countess™ Automated Cell Counter to determine the viability and total number of cells; the cells should have more than doubled in number.
- Carefully aspirate most of the supernatant, leaving approximately 100–200 µL behind. Remove the remaining medium with a 200-µL pipette.
- Resuspend the cell pellet in Resuspension Buffer T (included with Neon™ Transfection kits) at a final concentration of  $1.0 \times 10^7$ – $1.4 \times 10^7$  cells/mL.  
**Note:** Buffer T is required for blood cell transfections.
- Transfer the cells (10 µL per transfection reaction) to a sterile 1.5-mL microcentrifuge tube.
- Turn on the Neon™ unit and enter the following electroporation parameters in the Input window.

Pulse voltage	Pulse width	Pulse number	Cell density	Tip type
1650 V	10 ms	3	$1 \times 10^7$ cells/mL	10-µL

- Fill the Neon™ Tube with 3 mL Electrolytic Buffer (use Buffer E for the 10-µL Neon™ Tip).
- Insert the Neon™ Tube into the Neon™ Pipette Station until you hear a click.
- Transfer 1 µL of Epi5™ Reprogramming Vectors to the tube containing cells.
- Transfer 1 µL of Epi5™ p53 & EBNA Vectors to the tube containing cells and mix gently

12. Insert a 10- $\mu$ L Neon™ Tip into the Neon™ Pipette.
13. Press the push-button on the Neon™ Pipette to the first stop and immerse the Neon™ Tip into the cell-DNA mixture. Slowly release the push-button on the pipette to aspirate the cell-DNA mixture into the Neon™ Tip.
 

**Note:** Avoid air bubbles during pipetting to avoid arcing during electroporation. If you notice air bubbles in the tip, discard the sample and carefully aspirate fresh sample into the tip again without any air bubbles.
14. Insert the Neon™ Pipette with the sample vertically into the Neon™ Tube placed in the Neon™ Pipette Station until you hear a click.
15. Ensure that you have entered the appropriate electroporation parameters and press **Start** on the Neon™ touchscreen to deliver the electric pulse.
 

**Note:** After the electric pulse is delivered, the touchscreen displays “Complete” to indicate that electroporation is complete.
16. Remove the Neon™ Pipette from the Neon™ Pipette Station and immediately transfer the samples from the Neon™ Tip into one well of the pre-warmed 6-well plate with MEF feeders and containing 2 mL of complete StemPro™-34 medium with cytokines.
 

**Note:** Evenly distribute cells over the well in a dropwise manner.
17. Discard the Neon™ Tip into an appropriate biological hazardous waste container.
18. Repeat the process for any additional samples. Do not use Neon™ tip more than twice.
19. Incubate the plates at 37°C in a humidified CO<sub>2</sub> incubator overnight.

**Day 1:  
Switch to N2B27  
Medium**

1. Do not aspirate the spent medium from the plates. Add 1 mL of N2B27 Medium supplemented with 100 ng/mL bFGF (added freshly prior to use; see page 28 for recipe) to each well.

**Day 2 to 6:  
Add fresh N2B27  
Medium**

1. Carefully aspirate 0.5–1 mL of the supernatant from the top of the well, trying to not remove any of the suspension StemPro™ CD34<sup>+</sup> cells while they are still attaching.
2. Add 0.5–1 mL of N2B27 medium supplemented with 100 ng/mL bFGF.

**Day 7 to 9:  
Replace spent  
N2B27 Medium**

1. Carefully aspirate the spent medium and replace with 2 mL of N2B27 medium supplemented with 100 ng/mL bFGF.
2. Replace spent medium with 2 mL fresh medium every day, up to day 9 post-transfection.

**Day 9:  
Switch to human  
iPSC Medium**

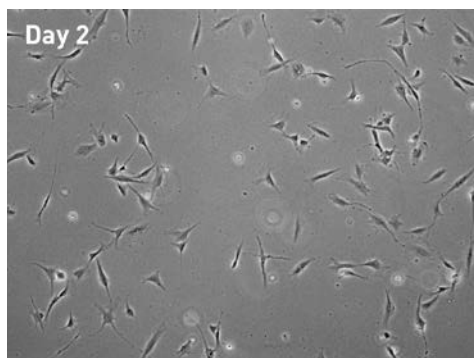
1. Aspirate the spent N2B27 medium and replace it with complete human iPSC Medium (see page 29 for recipe). Resume medium changes every day.
2. Observe the plates every other day under a microscope for the emergence of cell clumps indicative of transformed cells (see **Identify iPSC colonies**, page 19). Within 15 to 21 days of transfection, the iPSC colonies will grow to an appropriate size for transfer.

## Identify iPSC colonies

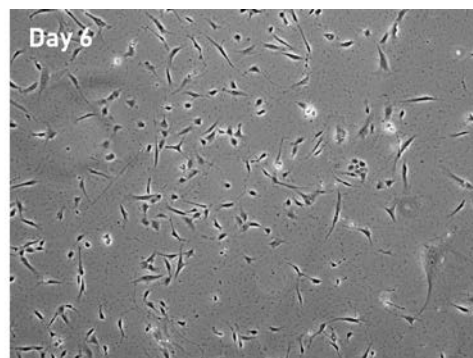
---

**Visual identification** By day 21, the colonies on the plates will appear large and compact and cover the majority of the surface area of the culture vessel. However, only a fraction of these colonies will consist of iPSCs, which exhibit a hESC-like morphology characterized by a flatter cobblestone-like appearance with individual cells clearly demarcated from each other in the colonies (see **Morphology of reprogrammed cells**, page 20). Therefore, we recommend that you perform live surface marker staining with Tra1-60 or Tra1-81 antibodies that recognize undifferentiated iPSCs (see **Live stain with antibodies** page 21). Note that other methods of identifying iPSCs, such as alkaline phosphatase staining, are also acceptable.

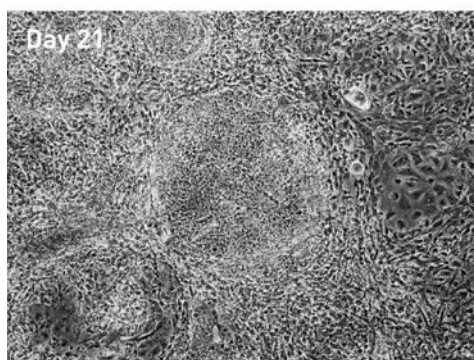
The images below illustrate the expected morphological changes as iPSC colonies emerge from human neonatal foreskin fibroblast cells (strain BJ) that were reprogrammed using the Epi5™ Episomal iPSC Reprogramming Kit. The phase contrast images were obtained at 5× magnification on Day 2, Day 6, and Day 21 post-transfection.



Day 2  
Visible morphology changes; cells appear unhealthy.



Day 6  
Groups of cube/round-shaped cells become visible throughout the plate

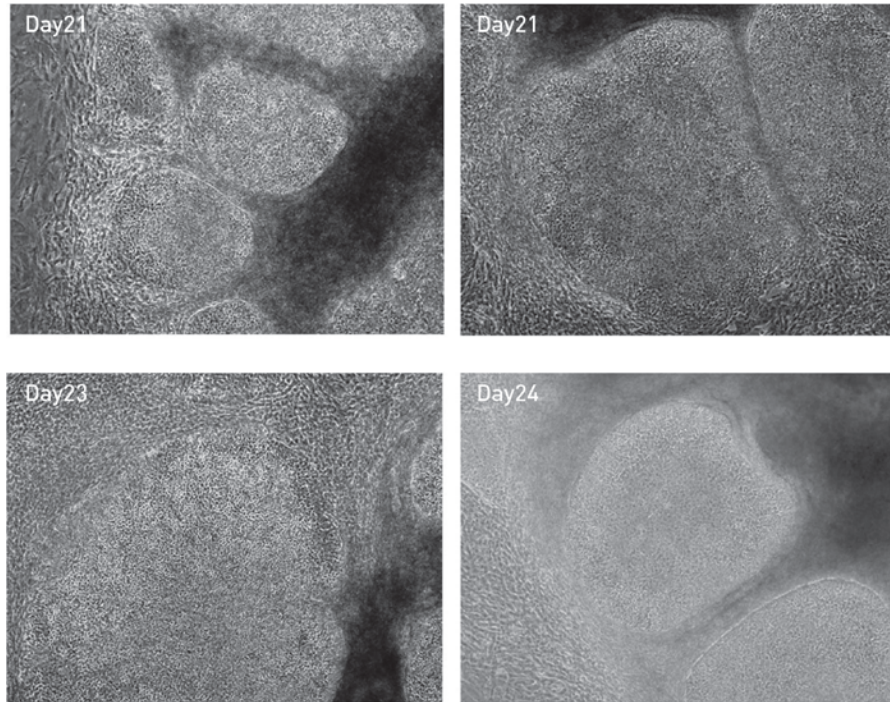


Day 21  
iPSC colonies emerge.

**Note:** Although colonies of “transformed” cells may emerge as early as 7 days after transfection, most of these colonies will not be correctly “reprogrammed” cells. iPSCs usually emerge a little later (around day 14 post-transfection), resemble embryonic stem cells in morphology, and express the cell surface markers Tra1-60 and Tra1-81.

## Morphology of reprogrammed cells

The images below show the expected morphology of emerging iPSCs generated by episomal reprogramming of human neonatal foreskin fibroblast cells (strain BJ). In these 5X images, lots of large, nested colonies are visible.



## Live stain with antibodies

---

One of the fastest and most reliable methods for selecting a reprogrammed colony is live staining with Tra1-60 or Tra1-81 antibodies that recognize undifferentiated iPSCs and enable the identification of reprogrammed cells from a variety of human cell types. Note that other methods of identifying iPSCs (such as alkaline phosphatase staining) are also acceptable.

**Required antibodies** Choose from the following:

- TRA-1-60 Alexa Fluor™ 488 Conjugate Kit for Live Cell Imaging (Cat. no. A25618)
- TRA-1-60 Alexa Fluor™ 594 Conjugate Kit for Live Cell Imaging (Cat. No. A24882)
- TRA-1-60 Alexa Fluor™ 555 Conjugate Kit for Live Cell Imaging (Cat. No. A24879)

**Live staining protocol**

**Caution:** Use aseptic technique when handling live-cell imaging reagents to prevent contamination.

1. Centrifuge the dye-conjugated antibody solution (e.g., 2 minutes at 10,000 × g) and only use the supernatant.  
**Note:** This step minimizes transferring protein aggregates that may have formed during storage, thereby reducing non-specific background staining.
2. Add a 1:50 volume of the dye-conjugated antibody directly to the cell culture medium of the cells to be stained (e.g., for each well of a 6-well plate, 10 uL of antibody and 1 mL of medium). Mix by gentle swirling.
3. Incubate for 30 minutes at 37°C.
4. Remove the staining solution and gently wash the cells 2–3 times with FluoroBrite™ DMEM.
5. For optimal results, image the cells immediately (i.e., within 30 minutes).  
**Note:** To continue culturing the cells, replace the FluoroBrite™ DMEM with fresh cell culture medium and return the cells to the 37°C incubator.

**Expected results**

Successful antibody staining can very specifically distinguish reprogrammed colonies from just plain transformed counterparts, and can be detected for up to 24–36 hours. This is particularly useful because it helps identifying and tracking of candidate iPS colonies before picking and the day after they are transferred into a new culture dish for expansion.

## Pick iPSC colonies

---

### Protocol for picking iPSC colonies (feeder-free culture)

1. Place the culture dish containing the reprogrammed cells under an inverted microscope and examine the colonies under 10X magnification.
2. Mark the colony to be picked on the bottom of the culture dish.  
**Note:** We recommend picking at least 10 distinct colonies by the end of each reprogramming experiment and expanding them in separate Geltrex™ matrix-coated 12-well culture plates.
3. Transfer the culture dish to a sterile cell culture hood (i.e., biosafety cabinet) equipped with a stereomicroscope.
4. Using a 25 gauge 1½ inch needle, cut the colony to be picked into 10–30 pieces in a grid-like pattern.
5. Using a 200-µL pipette, transfer the cut pieces to a freshly prepared Geltrex™ matrix-coated 12-well plate containing Essential 8™ Medium.
6. Incubate the Geltrex™ matrix-coated-coated plate containing the picked colonies in a 37°C, 5% CO<sub>2</sub> incubator.
7. Allow the colonies to attach to the culture plate for 48 hours before replacing the spent medium with fresh Essential 8™ Medium. After that, change the medium every day.
8. Treat the reprogrammed colonies like normal human iPSC colonies; passage, expand, and maintain them using standard culture procedures until you have frozen cells from two 60-mm plates (see page 33).

**Note:** Cells cultured in Essential 8™ Medium on Geltrex™ matrix-coated culture vessels should be passaged using 0.5 mM EDTA in DPBS. Use of enzymes such as collagenase and dispase for passaging these cells results in compromised viability and attachment.

**Note:** Newly derived iPSC lines may contain a fair amount of differentiation through passage 4. **It is not necessary to remove differentiated material prior to passaging.** By propagating/splitting the cells the overall culture health should improve throughout the early passages.

**Protocol for picking iPSC colonies (feeder-dependent)**

1. Place the culture dish containing the reprogrammed cells under an inverted microscope and examine the colonies under 10X magnification.
2. Mark the colony to be picked on the bottom of the culture dish.  
**Note:** We recommend picking at least 10 distinct colonies by the end of each reprogramming experiment and expanding them in separate 12-well MEF culture plates.
3. Transfer the culture dish to a sterile cell culture hood (i.e., biosafety cabinet) equipped with a stereomicroscope.
4. Using a 25 gauge 1½ inch needle, cut the colony to be picked into 10–30 pieces in a grid-like pattern.
5. Using a 200 µL pipette, transfer the cut pieces to a freshly prepared 12-well MEF culture plate (see page 31) containing human iPSC medium (see page 29).
6. Incubate the MEF culture plate containing the picked colonies in a 37°C incubator with a humidified atmosphere of 5% CO<sub>2</sub>.
7. Allow the colonies to attach to the culture plate for 48 hours before replacing the spent medium with fresh human iPSC medium. After that, change the medium every day.
8. Treat the reprogrammed colonies like normal human ESC colonies and passage, expand, and maintain them using standard culture procedures until you have frozen cells from two 60-mm plates (see page 34).

**Adapt iPSCs to feeder-free culture**

iPSCs cultured on MEF feeder layers can be adapted to feeder-free conditions in Essential 8™ Medium.

1. Harvest PSCs using an enzymatic method, such as Collagenase Type IV (Cat. no. 17104-019) or Dispase (Cat. no. 17105-041).
2. Plate cells on Geltrex™ Matrix-coated plates in Essential 8™ Medium (see page 30 for coating plates with Geltrex™ Matrix).
3. Incubate cells at 37°C, 5% CO<sub>2</sub> overnight.
4. Feed cells with Essential 8™ Medium beginning on the second day after splitting. Replace spent medium daily.

**Note:** Newly derived iPSC lines may contain a fair amount of differentiation through passage 4. **It is not necessary to remove differentiated material prior to passaging.** By propagating/splitting the cells the overall culture health should improve throughout the early passages.

## Detect episomal vectors by PCR

### Harvest iPSCs for PCR (feeder-free culture)

- Aspirate the medium from the dish containing iPSCs with a Pasteur pipette, and rinse the dish twice with Dulbecco's PBS (DPBS) without Calcium and Magnesium. Refer to table below for the recommended volumes.
- Add 0.5 mM EDTA in DPBS to the dish containing iPSCs. Adjust the volume of EDTA for various dish sizes. Swirl the dish to coat the entire cell surface.
- Incubate the vessel at room temperature for 5–8 minutes or 37°C for 4–5 minutes. When the cells start to separate and round up, and the colonies will appear to have holes in them when viewed under a microscope, they are ready to be removed from the vessel.  
**Note:** In larger vessels or with certain cell lines, this may take longer than 5 minutes.
- Aspirate the EDTA solution with a Pasteur pipette and add pre-warmed complete Essential 8™ Medium to the dish according to the table below.
- Remove the cells by gently squirting the colonies from the well using a 5-mL glass pipette. Avoid creating bubbles. Collect cells in a 15-mL conical tube.  
**Note: Do not scrape the cells from the dish.** There may be obvious patches of cells that were not dislodged and left behind. Do not attempt to recover them through scraping.  
**Note:** Depending upon the cell line, work with no more than one to three wells at a time, and work quickly to remove cells after adding Essential 8™ Medium to the well(s). The initial effect of the EDTA will be neutralized quickly by the medium. Some lines re-adhere very rapidly after medium addition, and must be removed 1 well at a time. Others are slower to re-attach, and may be removed 3 wells at a time.
- Centrifuge the cell suspension at 200 × g for 5 minutes to pellet cells.
- Aspirate and discard the supernatant. Resuspend the cell pellet in 500 µL DPBS and transfer resuspended cells to a thin-walled 0.5-mL PCR tube.
- Centrifuge the cell suspension at 200 × g for 5 minutes to pellet cells.
- Aspirate and discard the supernatant. Resuspend the cell pellet in 20 µL of Resuspension Buffer with 2 µL of Lysis Solution added to the Resuspension Buffer.
- Incubate the cells for 10 minutes in an incubator or thermal cycler that has been preheated to 75°C.
- Centrifuge the tube briefly to collect any condensation. Use 3 µL of the cell lysate in a 50-µL PCR reaction (see page 26).

Culture Vessel	Approximate Surface Area (cm <sup>2</sup> )	DPBS (mL)	0.5 mM EDTA in DPBS (mL)	Complete Essential 8™ Medium (mL)
6-well plate	10 cm <sup>2</sup> /well	2 mL/well	1 mL/well	2 mL/well
12-well plate	4 cm <sup>2</sup> /well	1 mL/well	0.4 mL/well	1 mL/well
24-well plate	2 cm <sup>2</sup> /well	0.5 mL/well	0.2 mL/well	0.5 mL/well
35-mm dish	10 cm <sup>2</sup>	2 mL	1 mL	2 mL
60-mm dish	20 cm <sup>2</sup>	4 mL	2 mL	4 mL
100-mm dish	60 cm <sup>2</sup>	12 mL	6 mL	12 mL



**Harvest iPSCs for PCR (feeder-dependent culture)**

1. Aspirate the MEF medium from a dish containing inactivated MEFs and add pre-warmed human iPSC Medium to the dish, 3–4 hours before harvesting hESCs.
2. Aspirate the spent medium from the dish with a Pasteur pipette.
3. Add Collagenase Type IV (1 mg/mL) solution to the dish containing hESCs. Adjust the volume of Collagenase Type IV for various dish sizes (e.g., 35-mm dishes require 1 mL of Collagenase IV).
4. Incubate the dish(es) for 30–60 minutes in a 37°C, 5% CO<sub>2</sub> incubator. Note that the incubation times may vary among different batches of collagenase; therefore, examination of the colonies is needed to determine the appropriate incubation time.  
**Note:** As an alternative to Collagenase Type IV, you may use 2 mg/mL of Dispase and incubate the dish for 2–3 minutes in a 37°C, 5% CO<sub>2</sub> incubator.
5. Stop the incubation when the edges of the colonies are starting to pull away from the plate.
6. Aspirate the Collagenase Type IV Solution with a Pasteur pipette. Remove the collagenase carefully without disturbing the attached cell layer.
7. Add human iPSC Medium to the dish. Use a 5-mL pipette to gently blow the cells off the surface of the dish while pipetting up and down. Make sure to pipet gently to minimize the formation of bubbles.
8. After the cells have been removed from the surface of the well, pool the contents of the wells into a 15-mL conical tube.
9. Using a 5-mL pipette, add human iPSC Medium to the dish to wash and collect any residual cells. Pipet up the medium and cells, and then add the collected cells to the 15-mL tube.
10. Pipet cells up and down gently a few times in the 15-mL tube to further break up cell colonies. Pipet carefully to reduce foaming.  
**Note:** Avoid making a single cell suspension.
11. Centrifuge the cell suspension at 200 × g for 5 minutes to pellet cells.
12. Aspirate and discard the supernatant. Resuspend the cell pellet in 500 µL DPBS and transfer resuspended cells to a thin-walled 0.5-mL PCR tube.
13. Centrifuge the cell suspension at 200 × g for 5 minutes to pellet cells.
14. Aspirate and discard the supernatant. Resuspend the cell pellet in 20 µL of Resuspension Buffer with 2 µL of Lysis Solution added to the Resuspension Buffer.
15. Incubate the cells for 10 minutes in an incubator or thermal cycler that has been preheated to 75°C.
16. Centrifuge the tube briefly to collect any condensation. Use 3 µL of the cell lysate in a 50-µL PCR reaction (see page 26).

**PCR primers for detecting Epi5™ Episomal Vectors**

You can detect the presence of Episomal iPSC Reprogramming Vectors in reprogrammed iPSC colonies by endpoint PCR using the PCR primers listed below. The EBNA-1 primer set can detect all five episomal plasmids in the kit. The oriP primer set can detect all episomal plasmids in the kit except pCXB-EBNA1, which lacks the OriP gene.

Transgene	Primers	Sequence	Expected size
oriP	pEP4-SF1-oriP	5'-TTC CAC GAG GGT AGT GAA CC-3'	544 bp
	pEP4-SR1-oriP	5'-TCG GGG GTG TTA GAG ACA AC-3'	
EBNA-1	pEP4-SF2-oriP	5'-ATC GTC AAA GCT GCA CAC AG-3'	666 bp
	pEP4-SR2-oriP	5'-CCC AGG AGT CCC AGT AGT CA-3'	

**PCR using AccuPrime™ High Fidelity Taq DNA Polymerase**

1. Add the components listed below to a DNase/RNase-free, thin-walled PCR tube. For multiple reactions, prepare a master mix of common components to minimize reagent loss and enable accurate pipetting.

**Note:** Assemble PCR reactions in a DNA-free environment. We recommend use of clean dedicated automatic pipettors and aerosol resistant barrier tips.

Component	Volume per reaction
10X PCR Buffer II	5 µL
Forward primer (10 µM stock)	1 µL
Reverse primer (10 µM stock)	1 µL
AccuPrime™ Taq Polymerase (5 units/µL)	1 µL
Cell Lysate	3 µL
Sterile distilled water	39 µL

2. Cap the tube, tap gently to mix, and centrifuge briefly to collect the contents.
3. Place the tube in the thermal cycler and use the PCR parameters shown below.

Step	Temperature	Time	Cycles
Initial denaturation	94°C	2 minutes	—
Denaturation	94°C	30 seconds	35–40
Annealing	55°C	30 seconds	
Elongation	72°C	1 minute	
Final elongation	72°C	7 minutes	—

4. Analyze the PCR products using 2% agarose gel electrophoresis.

# Appendix A: Recipes

## Media and reagents

### Basic FGF stock solution

- To prepare 10 mL of 10- $\mu$ g/mL Basic FGF solution, aseptically mix the following components:

Basic FGF	100 $\mu$ g
DPBS without Ca <sup>2+</sup> and Mg <sup>2+</sup>	9.8 mL
7.5% BSA	133 $\mu$ L
- Aliquot and store the Basic FGF solution at  $-20^{\circ}\text{C}$  for up to 6 months.  
**Note:** If desired, you may use 100  $\mu$ L of KnockOut™ Serum Replacement (KSR) instead of 7.5% BSA.

### 0.5 mM EDTA in DPBS

- To prepare 50 mL of 0.5 mM EDTA in DPBS, aseptically mix the following components in a 50-mL conical tube:

DPBS without Ca <sup>2+</sup> and Mg <sup>2+</sup>	50 mL
0.5 M EDTA	50 $\mu$ L
- Filter sterilize the solution and store at room temperature for up to 6 months.

### 1 mg/mL Collagenase Type IV Solution

- Add DMEM/F-12 to Collagenase Type IV to make a 10 mg/mL stock solution. Gently vortex to suspend, and filter sterilize the solution. This solution can be aliquoted and frozen at  $-20^{\circ}\text{C}$  until use.
- Make a working solution of 1 mg/mL Collagenase Type IV in DMEM/F-12. The working solution can be used for 2 weeks if properly stored at  $2-8^{\circ}\text{C}$  (store in aliquots to avoid repeated warming).

### Fibroblast Medium

To prepare 100 mL of complete Fibroblast Medium, aseptically mix the following components. Complete Fibroblast Medium can be stored at  $2-8^{\circ}\text{C}$  for up to 1 week.

Component	Volume
DMEM with GlutaMAX™-I	89 mL
FBS, ESC Qualified	10 mL
MEM Non-Essential Amino Acids Solution (10 mM)	1 mL

## N2B27 Medium

To prepare 250 mL of complete N2B27 Medium, aseptically mix the following components. N2B27 Medium (without bFGF) can be stored at 2–8°C for up to 1 week.

Component	Volume
DMEM/F-12 with HEPES	238.75 mL
N-2 Supplement (100X)	2.5 mL
B-27™ Supplement (50X)	5.0 mL
MEM Non-essential Amino Acids Solution (10 mM)	2.5 mL
GlutaMAX™-I (100X)	2.5 mL
β-Mercaptoethanol	454 µL
Basic FGF* (10 µg/mL)	2.5 mL

\* Prepare the N2B27 Medium without bFGF, and then supplement with fresh bFGF to a final concentration of 100 ng/mL when the medium is used.

## Essential 8™ Medium

1. Thaw the frozen Essential 8™ Supplement at 2–8°C overnight, or for 1–2 hours at room temperature. **Do not thaw the frozen supplement at 37°C.**
2. Mix the thawed supplement by gently inverting the vial a couple of times, remove 10 mL from the bottle of DMEM/F-12 (HAM) 1:1, and then aseptically transfer the entire contents of the Essential 8™ Supplement to the bottle of DMEM/F-12 (HAM) 1:1.
3. Swirl the bottle to mix and to obtain 500 mL of homogenous complete medium.
4. Complete Essential 8™ Medium can be stored at 2–8°C for up to 2 weeks. Before use, warm complete medium required for that day at room temperature until it is no longer cool to the touch. **Do not warm the medium at 37°C.**

## StemPro™-34 Medium

1. Thaw the frozen StemPro™-34 Nutrient Supplement at 2–8°C overnight.
2. After thawing, mix the supplement well by gently inverting the vial a couple of times, and then aseptically transfer the entire contents of the vial to the bottle of StemPro™-34 SFM. Swirl the bottle to mix and to obtain a homogenous medium.
3. Aseptically add GlutaMAX™-I (100X) supplement (Cat. no. 25030) to a final concentration of 1X (5 mL of 100X GlutaMAX™-I supplement to 500 mL of medium).
4. Before use, add the following cytokines to the indicated final concentration into the medium. **These must be added freshly, just prior to use.**
  - Recombinant human SCF 100 ng/mL
  - Recombinant human IL-3 50 ng/mL
  - Recombinant human GM-CSF 25 ng/mL
5. Complete StemPro™-34 Medium (without the cytokines) can be stored at 2–8°C **in the dark** for up to 4 weeks. The thawed nutrient supplement has a shelf life of 2 weeks when stored at 2–8°C, in the dark.

## Human iPSC Medium

To prepare 100 mL of Human iPSC Medium, aseptically mix the components listed below. Human iPSC Medium can be stored at 2–8°C for up to 1 week.

Component	Volume
DMEM/F-12 (1X) with GlutaMAX™-I	79 mL
MEM Non-essential Amino Acids Solution (10 mM)	1 mL
β-Mercaptoethanol (1000X)	100 μL
KnockOut™ Serum Replacement (KSR)	20 mL
Basic FGF* (10 μg/mL)	1 mL

\* Prepare the iPSC Medium without bFGF, and then supplement with fresh bFGF to a final concentration of 100 ng/mL when the medium is used.

## MEF Medium

To prepare 100 mL of complete MEF medium, aseptically mix the components listed below. Complete MEF Medium can be stored at 2–8°C for up to 1 week.

Component	Volume
DMEM with GlutaMAX™-I	89 mL
FBS, ESC-Qualified	10 mL
MEM Non-essential Amino Acids Solution (10 mM)	1 mL
β-Mercaptoethanol (1000X)	100 μL

# Appendix B: Prepare culture vessels

## Coat culture vessels with Geltrex™ matrix

### Coating protocol

1. Thaw a 5-mL bottle of Geltrex™ LDEV-Free hESC-Qualified Reduced Growth Factor Basement Membrane Matrix at 2–8°C overnight.
2. For reprogramming experiments using fibroblasts, dilute the thawed Geltrex™ matrix solution 1:1 with cold sterile DMEM/F-12 to prepare 1-mL aliquots (or another volume suitable for your needs) in tubes chilled on ice. These aliquots can be frozen at –20°C or used immediately.  
  
For reprogramming experiments using hemopoietic progenitor cells (HPCs), such as StemPro™ CD34<sup>+</sup> cells, do **not** dilute the Geltrex™ matrix solution 1:1 to create the intermediate dilution. The Geltrex™ matrix solution needs to be twice as strong for reprogramming experiments with HPCs.
3. To create working stocks, dilute an aliquot of Geltrex™ matrix solution 1:50 with cold DMEM on ice. This creates a final dilution of 1:100 for fibroblasts or a final dilution of 1:50 for StemPro™ CD34<sup>+</sup> cells.  
**Note:** An optimal dilution of the Geltrex™ matrix solution may need to be determined for each cell line. Try various dilutions from 1:30 to 1:100.
4. Quickly cover the whole surface of each culture dish with the Geltrex™ matrix solution (see table below).
5. Incubate the dishes in a 37°C, 5% CO<sub>2</sub> incubator for 1 hour.
6. Geltrex™ matrix-coated culture dishes can now be used or stored at 2–8°C for up to a week. Do not allow dishes to dry.
7. Aspirate the diluted Geltrex™ matrix solution from the culture dish and discard. You do not need to rinse off the Geltrex™ matrix solution from the culture dish after removal. Cells can now be passaged directly onto the Geltrex™ matrix-coated culture dish.

Culture vessel	Surface area	Volume of Geltrex™ matrix dilution
6-well plate	10 cm <sup>2</sup> /well	1.5 mL/well
12-well plate	4 cm <sup>2</sup> /well	750 µL/well
24-well plate	2 cm <sup>2</sup> /well	350 µL/well
35-mm dish	10 cm <sup>2</sup>	1.5 mL
60-mm dish	20 cm <sup>2</sup>	3.0 mL
100-mm dish	60 cm <sup>2</sup>	6.0 mL

## MEF culture dishes

---

### Gelatin coat culture vessels

1. Cover the whole surface of each culture vessel with Attachment Factor (AF) solution and incubate the vessels for 30 minutes at 37°C or for 2 hours at room temperature.  
**Note:** AF (Cat. no. S-006-100) is a sterile 1X solution containing 0.1 % gelatin available from Life Technologies (see page 36 for ordering information).
2. Using sterile technique in a laminar flow culture hood, completely remove the AF solution from the culture vessel by aspiration.  
**Note:** It is not necessary to wash the culture surface before adding cells or medium. Coated vessels may be used immediately or wrapped in Parafilm™ sealing film and stored at room temperature for up to 24 hours.

### Thaw MEFs

1. Remove the cryovial containing inactivated MEFs from the liquid nitrogen storage tank.
2. Briefly roll the vial between hands to remove frost, and swirl it gently in a 37°C water bath.
3. When only a small ice crystal remains in the vial, remove it from water bath. Spray the outside of the vial with 70% ethanol before placing it in the cell culture hood.
4. Pipet the thawed cells gently into a 15-mL conical tube.
5. Rinse the cryovial with 1 mL of pre-warmed MEF medium (see page 29). Transfer the medium to the same 15-mL tube containing the cells.
6. Add 4 mL of pre-warmed MEF medium **dropwise** to the cells. Gently mix by pipetting up and down.  
**Note:** Adding the medium slowly helps the cells to avoid osmotic shock.
7. Centrifuge the cells at  $200 \times g$  for 5 minutes.
8. Aspirate the supernatant and resuspend the cell pellet in 5 mL of pre-warmed MEF medium.
9. Remove 20  $\mu$ L of the cell suspension and determine the viable cell count using your method of choice (e.g., Countess™ Automated Cell Counter).

### Plate MEFs

1. Centrifuge the remaining cell suspension (step 9, previous page) at  $200 \times g$  for 5 minutes at room temperature.
2. Aspirate the supernatant. Resuspend the cell pellet in MEF medium (see page 29) to a density of  $2.5 \times 10^6$  cells/mL.
3. Aspirate the gelatin solution from the gelatin coated culture vessel.
4. Add the appropriate amount of MEF medium into each culture vessel (refer to the table below).
5. Into each of these culture vessels, add the appropriate amount of MEF suspension (refer to the table below).  
**Note:** The recommended plating density for Gibco™ Mouse Embryonic Fibroblasts (Irradiated) (Cat. no. S1520-100) is  $2.5 \times 10^4$  cells/cm<sup>2</sup>.
6. Move the culture vessels in several quick back-and-forth and side-to-side motions to disperse the cells across the surface of the vessels.
7. Incubate the cells in a 37°C incubator with a humidified atmosphere of 5% CO<sub>2</sub>.
8. Use the MEF culture vessels within 3–4 days after plating.

Vessel size	Approximate growth area	Volume of MEF medium	Number of MEFs	Volume of MEF suspension
24-well plate	2 cm <sup>2</sup> /well	0.5 mL	5.0 × 10 <sup>4</sup> /well	20 μL
12-well plate	4 cm <sup>2</sup> /well	1 mL	1.0 × 10 <sup>5</sup> /well	40 μL
6-well plate	10 cm <sup>2</sup> /well	2 mL	2.5 × 10 <sup>5</sup> /well	0.1 mL
60-mm dish	20 cm <sup>2</sup>	5 mL	5.0 × 10 <sup>5</sup>	0.2 mL
100-mm dish	60 cm <sup>2</sup>	10 mL	1.5 × 10 <sup>6</sup>	0.6 mL
25-cm <sup>2</sup> flask	25 cm <sup>2</sup>	5 mL	6.3 × 10 <sup>5</sup>	0.25 mL
75-cm <sup>2</sup> flask	75 cm <sup>2</sup>	15 mL	1.9 × 10 <sup>6</sup>	0.75 mL



# Appendix C: Cryopreserve iPSCs

## Freeze iPSCs (Feeder-Free Culture)

### Freeze feeder-free iPSC cultures

1. Pre-warm the required volume of Essential 8™ Medium at room temperature until it is no longer cool to the touch. **Do not warm the medium in a 37°C water bath.**
2. Prepare Essential 8™ Freezing Medium. For every 1 mL of freezing medium needed, aseptically combine the following components in a sterile 15-mL tube:

Complete Essential 8™ Medium	0.9 mL
DMSO	0.1 mL
3. Place the tube with Essential 8™ Freezing Medium on ice until use. Discard any remaining freezing medium after use.
4. Aspirate the spent medium from the dish, and rinse the cells twice with DPBS without Ca<sup>2+</sup> and Mg<sup>2+</sup> (see table below).
5. Add 0.5 mM EDTA solution (see page 27) to the dish. Adjust the volume of EDTA for various dish sizes (see table below). Swirl the dish to coat the entire cell surface.
6. Incubate the dish at room temperature for 5–8 minutes or at 37°C for 4–5 minutes. When the cells start to separate and round up, and the colonies appear to have holes in them when viewed under a microscope, they are ready to be removed from the vessel.
7. Aspirate the EDTA solution with a Pasteur pipette.
8. Add ice-cold Essential 8™ Freezing Medium to the dish (see table below).
9. Remove the cells by gently squirting the colonies from the well using a 5-mL glass pipette. Avoid creating bubbles. Collect the cells in a 15-mL conical tube and place on ice.
10. Resuspend the cells gently. Aliquot 1 mL of the cell suspension into each cryovial.
11. Quickly place the cryovials in a cryofreezing container (e.g., Mr. Frosty) and freeze the cells by decreasing the temperature by 1°C per minute. Once frozen, transfer the cells to –80°C overnight.
12. After overnight storage at –80°C, transfer the cells to a liquid nitrogen tank vapor phase for long-term storage.

Culture vessel	Approximate surface area	DPBS (mL)	0.5 mM EDTA in DPBS (mL)	Essential 8™ Freezing Medium (mL)
6-well plate	10 cm <sup>2</sup> /well	2 mL/well	1 mL/well	1 mL/well
35-mm dish	10 cm <sup>2</sup>	2 mL	1 mL	1 mL
60-mm dish	20 cm <sup>2</sup>	4 mL	2 mL	2 mL
100-mm dish	60 cm <sup>2</sup>	12 mL	6 mL	6 mL

## Freeze iPSCs (Feeder-Dependent Culture)

---

### iPSC freezing medium for feeder-dependent culture

1. Prepare the Freezing Media A and B immediately before use.
2. In a sterile 15-mL tube, mix together the following reagents for every 1 mL of **freezing medium A** needed:

Human iPSC medium	0.5 mL
KnockOut™ Serum Replacement	0.5 mL
3. In another sterile 15-mL tube, mix together the following reagents for every 1 mL of **freezing medium B** needed:

Human iPSC medium	0.8 mL
DMSO	0.2 mL
4. Place the tube with freezing medium B on ice until use (you can keep freezing medium A at room temperature). Discard any remaining freezing medium after use.

### Freezing feeder-dependent iPSC cultures

1. Prepare the required volume of fresh freezing medium and place it on ice.
2. Aspirate the culture medium and rinse the dishes twice with DPBS without  $\text{Ca}^{2+}$  and  $\text{Mg}^{2+}$  (see table on page 35).
3. Gently add Collagenase IV solution (see page 27) to the culture dish. Adjust the volume of Collagenase IV solution for various dish sizes (see table on page 35).
4. Incubate the dish with cells for 5–20 minutes in a 37°C incubator with a humidified atmosphere of 5%  $\text{CO}_2$ .

**Note:** Incubation times may vary among different batches of collagenase. Therefore, the appropriate incubation time should be optimized by examining the colonies periodically under microscope during incubation.
5. Stop the incubation when the edges of the colonies are starting to pull away from the plate.
6. Remove the culture dish from the incubator, aspirate the Collagenase IV solution, and gently rinse the dish with DPBS without  $\text{Ca}^{2+}$  and  $\text{Mg}^{2+}$ .
7. Add human iPSC Medium or DMEM/F-12 (see table on page 35) and gently dislodge the cells off the surface of the culture dish using a sterile pipette or a cell scraper. Transfer the cells to a sterile 15-mL centrifuge tube. Rinse the dish with additional iPSC Medium or DMEM/F-12 to collect any leftover colonies.
8. Centrifuge the cells at  $200 \times g$  for 2–4 minutes at room temperature.
9. Discard the supernatant, gently tap the tube to dislodge the cell pellet from the tube bottom, and resuspend the cells in **freezing medium A** (see table on page 35). After the cell clumps have been uniformly suspended, add an equal volume of **freezing medium B** to the cell suspension in a drop-wise manner while gently swirling the cell suspension to mix.

**Note:** At this point, the cells are in contact with DMSO, and work must be performed efficiently with no or minimum delays. After the cells come into contact with DMSO, they should be aliquoted and frozen within 2–3 minutes.

10. Aliquot 1 mL of the cell suspension into each cryovial.
11. Quickly place the cryovials containing the cells in a cryo freezing container (e.g., Mr. Frosty) to freeze the cells at 1°C per minute and transfer them to -80°C overnight.
12. After overnight storage at -80°C, transfer the cells to a liquid nitrogen tank vapor phase for long term storage.

Culture Vessel	Approximate surface area	DPBS (mL)	Collagenase IV solution (mL)	Human iPSC medium (mL)	Freezing Medium A or B (mL)
6-well plate	10 cm <sup>2</sup> /well	2 mL/well	1 mL/well	1 mL/well	0.5 mL/well
35-mm dish	10 cm <sup>2</sup>	2 mL	1 mL	1 mL	0.5 mL
60-mm dish	20 cm <sup>2</sup>	4 mL	2 mL	2 mL	1 mL
100-mm dish	60 cm <sup>2</sup>	12 mL	6 mL	6 mL	3 mL

# Appendix D: Ordering information

## Accessory products

### Epi5™ episomal reprogramming products

For more information about the following products, refer to our website at [thermofisher.com](http://thermofisher.com) or contact Technical Support (see page 39).

Product	Quantity	Catalog No.
Epi5™ Episomal iPSC Reprogramming Kit	1 kit	A15960

### Media, sera, and supplements

For more information about the following products, refer to our website at [thermofisher.com](http://thermofisher.com) or contact Technical Support (see page 39).

Product	Quantity	Catalog No.
Dulbecco's Modified Eagle Medium (DMEM) with GlutaMAX™-I (high glucose)	500 mL	10569-010
DMEM/F-12 with HEPES	500 mL	11330-032
KnockOut™ DMEM/F-12	500 mL	12660-012
KnockOut™ Serum Replacement	500 mL	10828-028
Essential 8™ Medium	500 mL	A1517001
StemPro™-34 SFM (1X), Liquid	500 mL	10639-011
MEM Non-Essential Amino Acids Solution (10 mM)	100 mL	11140-050
Basic Fibroblast Growth Factor (bFGF), recombinant human	10 µg	PHG0264
Bovine Albumin Fraction V Solution (BSA), 7.5%	100 mL	15260-037
Fetal Bovine Serum (FBS), ES-Cell Qualified	500 mL	16141-079
β-Mercaptoethanol (1000X), liquid	50 mL	21985-023
GlutaMAX™-I Supplement	100 mL	35050-061
N-2 Supplement (100X)	5 mL	17502-048
B-27™ Serum-Free Supplement (50X)	10 mL	17504-044
DPBS without Ca <sup>2+</sup> or Mg <sup>2+</sup>	500 mL	14190-144
DMSO, anhydrous	10 × 3 mL	D12345

## Matrices and dissociation reagents

For more information about the following products, refer to our website at [thermofisher.com](http://thermofisher.com) or contact Technical Support (see page 39).

Product	Quantity	Catalog No.
Geltrex™ LDEV-Free hESC-Qualified Reduced Growth Factor Basement Membrane Matrix	5 mL	A1413302
Attachment Factor	100 mL	S-006-100
UltraPure™ 0.5 M EDTA, pH 8.0	4 × 100 mL	15575-020
0.05% Trypsin-EDTA (1X), Phenol Red	100 mL	25300-054
Collagenase, Type IV, powder	1 g	17104-019
Dispase II, powder	5 g	17105-041

## Cells

For more information about the following products, refer to our website at [thermofisher.com](http://thermofisher.com) or contact Technical Support (see page 39).

Product	Quantity	Catalog No.
Gibco™ Mouse Embryonic Fibroblasts (Irradiated)	1 mL	A24903
StemPro™ CD34+ Cell Kit (contains 0.5 × 10 <sup>6</sup> cells)	1 kit	A14059
Human Dermal Fibroblasts, neonatal (HDFn)	1 vial	C-004-5C
Human Dermal Fibroblasts, adult (HDFa)	1 vial	C-013-5C

## Antibodies

For more information about the following products, refer to our website at [thermofisher.com](http://thermofisher.com) or contact Technical Support (see page 39).

Product	Quantity	Catalog No.
TRA-1-60 Alexa Fluor™ 488 Conjugate Kit for Live Cell Imaging	50 tests	A25618
TRA-1-60 Alexa Fluor™ 594 Conjugate Kit for Live Cell Imaging	50 tests	A24882
TRA-1-60 Alexa Fluor™ 555 Conjugate Kit for Live Cell Imaging	50 tests	A24879

## Reagents for PCR detection of episomal vectors

For more information about the following products, refer to our website at [thermofisher.com](http://thermofisher.com) or contact Technical Support (see page 39).

Product	Quantity	Catalog No.
CellsDirect Resuspension & Lysis Buffers	1 kit	11739-010
AccuPrime™ Taq DNA Polymerase High Fidelity	200 reactions	12346-086

## Equipment

For more information about the following products, refer to our website at [thermofisher.com](http://thermofisher.com) or contact Technical Support (see page 39).

Product	Quantity	Catalog No.
Countess™ II Automated Cell Counter	1 unit	AMQAX1000
StemPro™ EZPassage™ Disposable Stem Cell Passaging Tool	10 units	23181-010
Neon™ Transfection System	1 unit	MPK5000
Neon™ Transfection System 10 µL Tip Kit	50 reactions	MPK1025
Neon™ Transfection System 100 µL Tip Kit	50 reactions	MPK10025

# Documentation and support

## Customer and technical support

---

Visit [www.thermofisher.com/support](http://www.thermofisher.com/support) for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
  - Product FAQs
  - Software, patches, and updates
- Order and web support
- Product documentation, including:
  - User guides, manuals, and protocols
  - Certificates of Analysis
  - Safety Data Sheets (SDS; also known as MSDSs)

**Note:** For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

### Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.lifetechnologies.com/termsandconditions](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

For support visit [thermofisher.com/support](http://thermofisher.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

[thermofisher.com](http://thermofisher.com)

07 March 2016

**ThermoFisher**  
S C I E N T I F I C