

# Hypoxia and iPSC: A Low Oxygen Atmosphere Supports the Xeno-free Generation, Expansion, and Differentiation of Human Induced Pluripotent Stem Cells

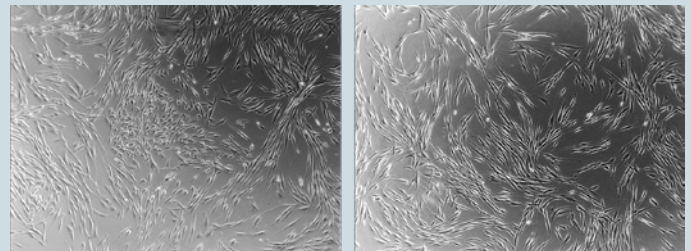
**Rick Cohen**  
Rutgers University, Piscataway, NJ, USA

## Abstract

We previously demonstrated that using low oxygen tension (4 %) increased the efficiency of reprogramming human somatic cells to pluripotency (Figure 1). In this study, we extend our findings and further increased the development of this culture paradigm using different coating substrates. We are able to observe that post-electroporated fibroblasts cultivated at 4 % O<sub>2</sub> grow on synthetic and biological surfaces. Further, we observed normal spreading of iPSC and a high level of purity in early iPSC passages. The cells were successfully differentiated into motor neuron and cardiomyocyte lineages emphasizing the efficacy of low oxygen levels during cell cultivation.

4 % O<sub>2</sub>

17 % O<sub>2</sub>



**Figure 1:** Generation of iPSCs under varying oxygen concentrations. After 12 days, early colony formation is seen in low O<sub>2</sub> cultures (left), whereas this is not observed as early in norm-oxy cultures (right).

[Read more in Application Note No. 338: Low Oxygen Levels Enhance the Efficiency of Reprogramming Human Somatic Cells to Pluripotency](#)

## Introduction

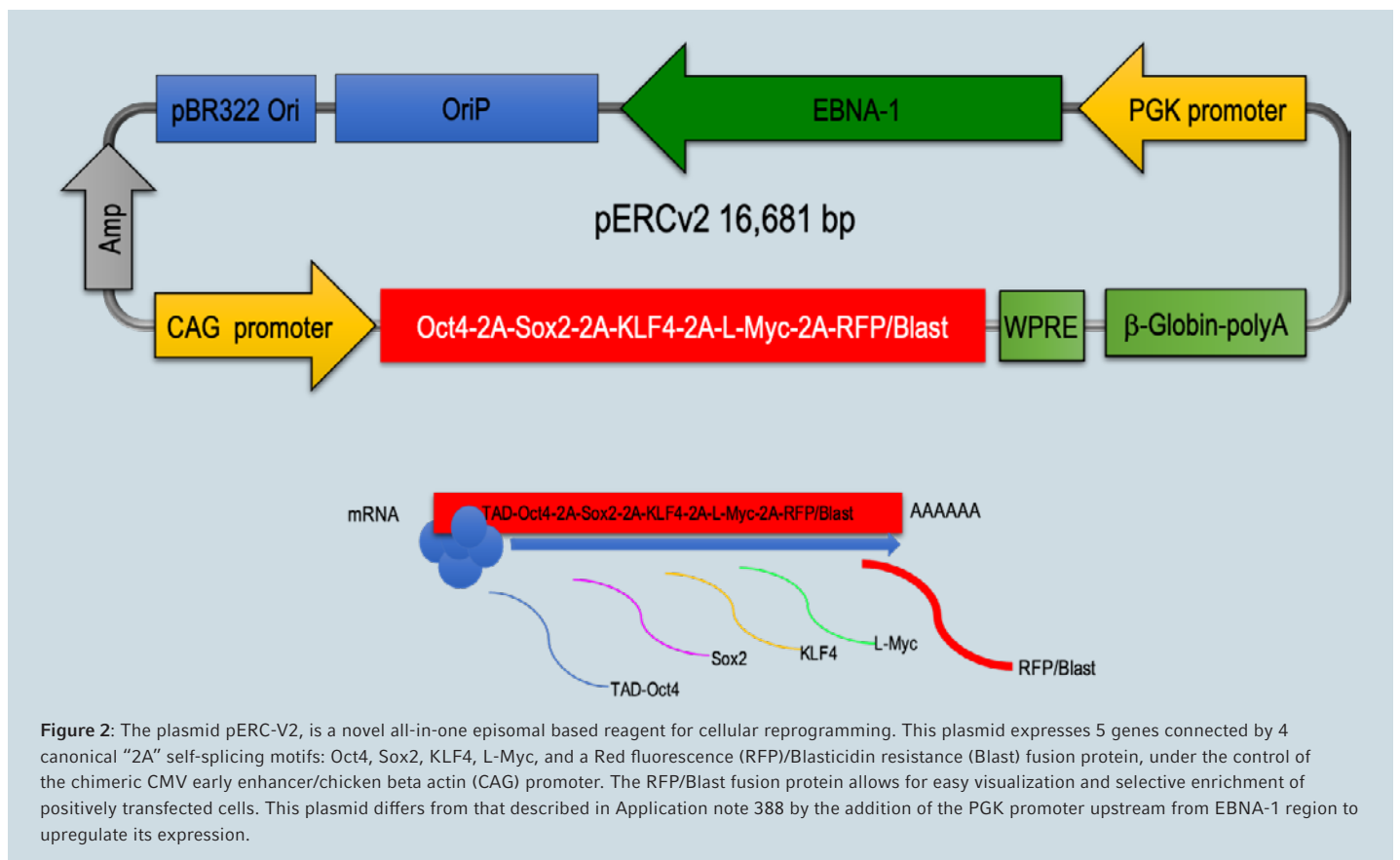
Reprogramming human somatic cells into the pluripotent state is a subject of thousands of publications since their discovery in 2007 [1, 2]. Originally these studies were pioneered with a set of four genes found to be expressed in native human embryonic stem cells; Oct4, Sox2, KLF4, and c-Myc (Lin28, and Nanog), delivered using genetically modifying methods such as retrovirus, and cultured on non-defined matrices. In the 13 years since these seminal studies many improvements were made such as (1) replacing genetically modifying methods with non-genome altering safer alternatives; (2) replacement of c-Myc oncogene with non-transforming family member, L-Myc; (3) inclusion of small molecules to boost efficiency of reprogramming; (4) optimizing culture conditions which

includes the use of low O<sub>2</sub> tension (4-5 %) and; (5) use of clinically relevant defined medias and matrices [3-11]. In this novel study, we combined many of the improvements to demonstrate the successful reprogramming of human foreskin fibroblasts with low O<sub>2</sub> conditions in the CellXpert® C170i CO<sub>2</sub> Incubator and an improved upon single fluorescent and selectable polycistronic Episomal vector as described in Application Note 338. We observed that the tested growth substrates offered stable cell adhesion and spreading following electroporation of fibroblasts, which in turn led to a robust production of karyotypically normal iPSC colonies capable of robust expansion in a defined media, and ability to differentiate into neural and cardiac lineages.

## Materials and Methods

The reagents and procedures used in this Application Note are similar to that described in Application Note 338. Notable differences or updates are summarized below, and more in-depth details are seen in Appendix 1. We have improved the pERCv1 plasmid as previously described by including a PGK promoter upstream from the EBNA-1 gene cassette (pERCv2, Figure 2, [12]). When indicated, culture vessels were coated with 5 ug/ml of Vitronectin as per manufacturer instructions (PeproTech®) and used in comparison with a synthetic substrate (Eppendorf FN1 motifs CCCadvanced® cultureware).

Once iPSC colonies appeared in Reprogramming Media, they were transitioned into Animal Free Low Protein hESC media for expansion (PeproTech, alternatively prepare as detailed in Chen et al., 2011 [13] or in Kuo et al., 2020 [14]). All iPSC expansion and early steps in neuronal differentiation was carried out with 6 well format dishes, whereas cells analyzed for immunostaining or terminally differentiated to motor neurons or cardiomyocytes were spin seeded (300 x g, 5 min at RT), onto 24 well format culture plates. Neurons were developed as described by Bianchi et al. 2018 [15] using the same base medias as described by Francis et al. 2019 [16]. Cardiomyocytes were generated using protocols as described by Burridge and Coworkers [17, 18].

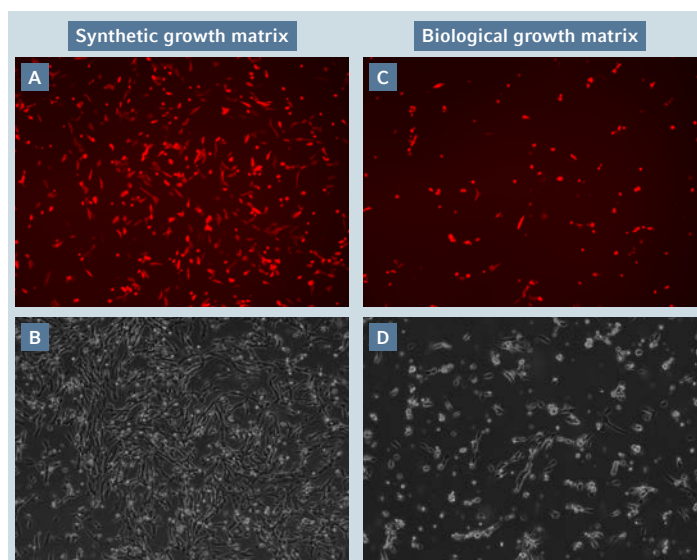


## Results and Discussion

### Reprogramming of Fibroblasts

Twenty four hours after electroporation, the fibroblast cultures plated onto biological and synthetic coatings were observed for red fluorescence (Figure 3). Even though a smaller number of cells ( $0.75 \times 10^6$  vs.  $1 \times 10^6$ ) were plated on the synthetic coating, it was clear that the cultures survived, and spread more robustly. This led to a failure to observe reprogramming in many of the Vitronectin plated cultures. We were able to observe reprogramming events as early as

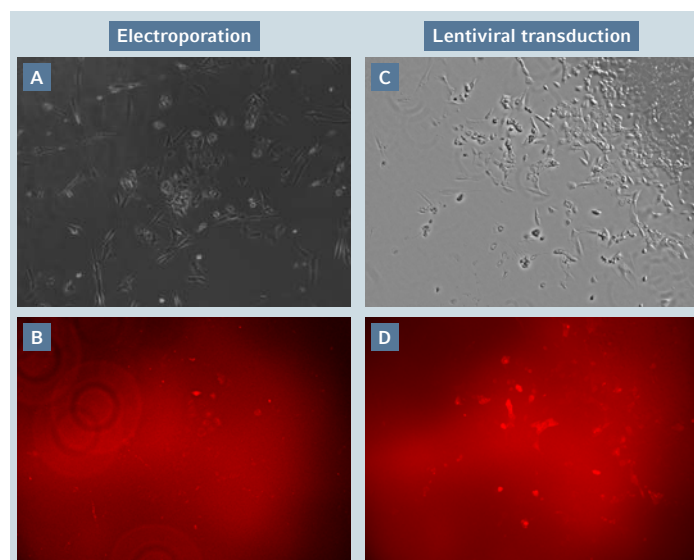
11 days post modification (Figure 4). The normally evenly distributed spindle shaped fibroblasts adopted a cuboidal appearance and organized into rudimentary colonies in both electroporated (Fig 4 A/B) and lentiviral transduced (Fig. 4 c-d) cultures. After 11 days only a minor amount of RFP was observed in both types of reprogramming paradigms, indicating possible loss of plasmid (episomal), or epigenetic inactivation (lentiviral).



**Figure 3:** RFP-expression 24h post electroporation. Fibroblasts were removed from 10 cm uncoated TC grade dishes using TrypLE, electroporated, and plated onto 6-well plates coated with synthetic ( $0.75 \times 10^6$  cells in A/B), or biological ( $1 \times 10^6$  cells in C/D) substrates.

### Expansion and Characterization of iPSCs under hypoxic conditions

The rudimentary colonies shown in figure 4 gave rise to more mature colonies at day 21 to day 30 post modification and were passaged using gentle non-enzymatic methods (Figure 5). After 3 routine passages, we observed that the colonies derived and cultured on the synthetic matrix

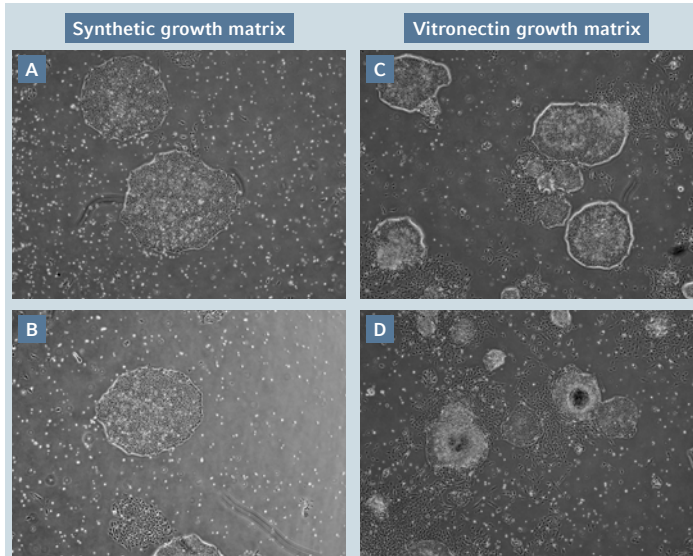


**Figure 4:** Appearance of presumptive reprogrammed colonies 11 days post electroporation (A/B) or lentiviral transduction (C/D). Fibroblasts ( $1 \times 10^6$ ) were electroporated as described or transduced with lentivirus containing the identical poly-cistronic supergene as seen in the Episomal plasmid. Cells under both conditions were cultured in a similar fashion. While RFP signal was significantly downregulated in both conditions (B/D), the 4 factors presumably reached a threshold for successful reprogramming at least by day 11 as seen in the appearance of small clusters (A/C). These small clusters went on to form mature looking colonies within 3 weeks.

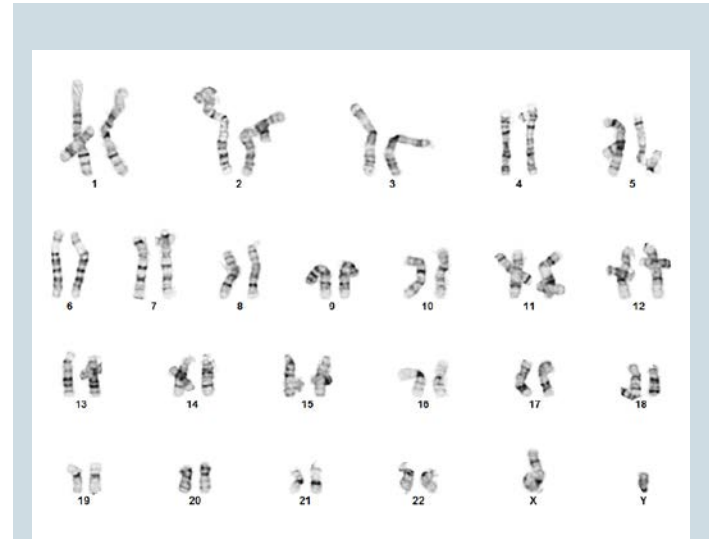
(A/B vs. C/D) appeared purer than on Vitronectin, with less appearance of non-reprogrammed fibroblasts. In order to purify the culture of Vitronectin derived cells, manual picking of colonies was used, whereas iPSCs grown on synthetic substrates outgrew any minor number of fibroblasts that were dislodged during the non-enzymatic passaging.

After 7 passages, the Episomal plasmid reprogrammed line from the synthetic coating was analyzed for karyotype and found to be normal (Figure 6). Similar cultures were spin seeded onto 24 well dishes and analyzed for routine pluripotency or differentiation markers. (Figure 7). At passage 8, the cultures were found to lack any appreciable expression

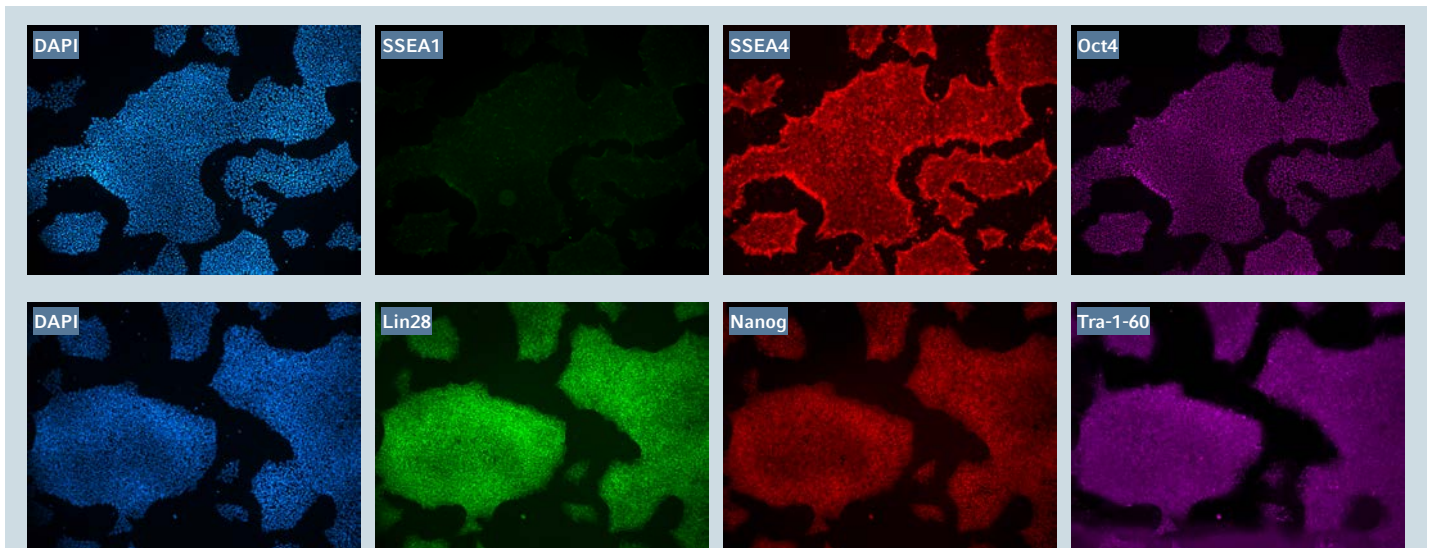
of SSEA1, a marker of differentiation, and robustly expressed both the SSEA4 surface marker, and Oct4 transcription factor. Likewise, these cell co-expressed Lin28 cytoplasmic RNA binding protein, Nanog transcription factor and the Tra-1-60 cell surface marker.



**Figure 5:** Standard passing techniques with iPSCs grown on different substrates. iPSCs grown on synthetic (A/B) or Vitronectin (C/D) matrices were developed and passaged using described techniques. After 3 routine passages using PBS-EGTA, the colonies developed on Vitronectin appeared to differentiate around the edges more readily and carry more non-reprogrammed or spontaneously differentiating fibroblasts into the serial cultures.



**Figure 6:** Karyotype analysis of PM1Ep5-ERCv2p7 iPSC Cells. Cytogenetic analysis was performed on 20 G-banded metaphase cells and demonstrated normal male karyotype.

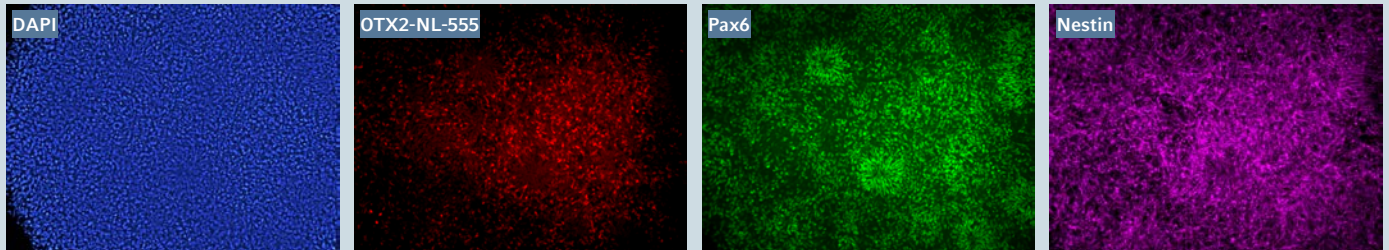


**Figure 7:** Immunofluorescence staining of markers for pluripotency and differentiation in low passage iPSC cells

### Differentiation of iPSCs to neuronal and cardiomyocyte lineages

Using standard published protocols, the cells were first differentiated into neuroepithelial cells (not shown) and then secondly into neural stem cells (Figure 8). After 4 passages

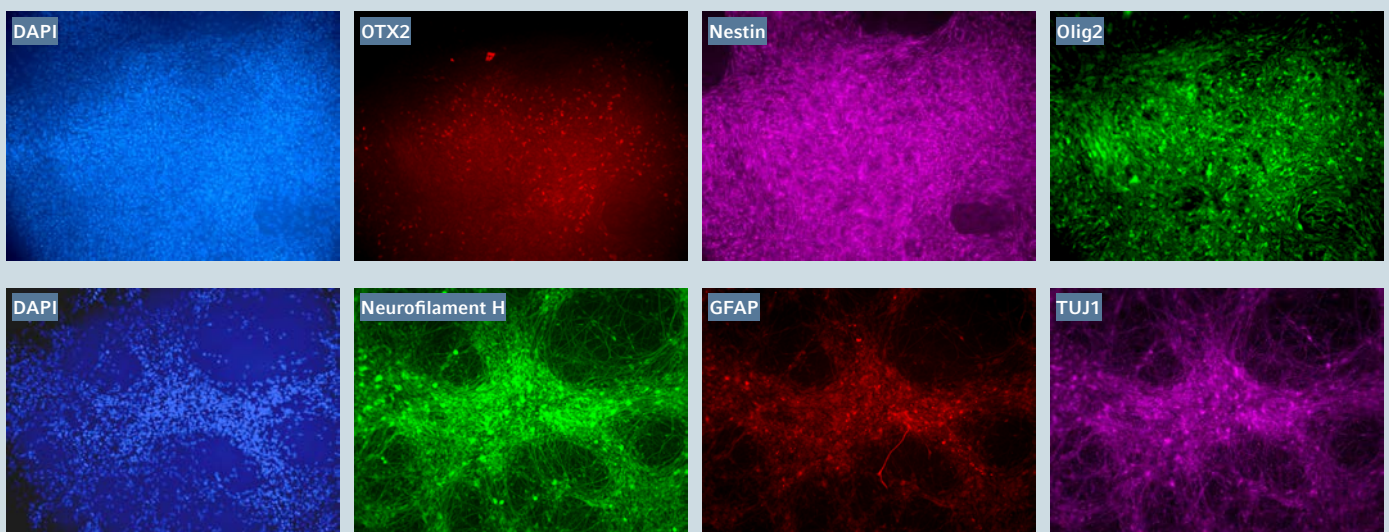
in neural stem cell media, the cells appeared to differentiate into a uniform mat (8A), with repeating patterns of rosettes (8C), characteristic of neural stem cell stage. Many of these cells were OTX2+, with a greater number expressing both Pax6, and Nestin.



**Figure 8:** Differentiation of iPSC into neural stem cells in low oxygen environment; iPSCs were differentiated using protocol as described by Bianchi and coworkers. Cells at second stage (MN2) were expanded using FGF-2 and EGF and are described as neural stem cells (NSC).

Since this was the second step out of four to derived motor neurons, we continued to differentiate the stem cells into presumptive motor neuron precursors. The small molecules and growth factors were altered to continue the differentiation protocol (Figure 9, upper panel), and some of the cells continued to express OTX2 and Nestin, while many began to express Olig2, a marker of motor neuron lineage. After one week under these conditions, the cells were finally exposed to a set of cytokines and small molecules that induced

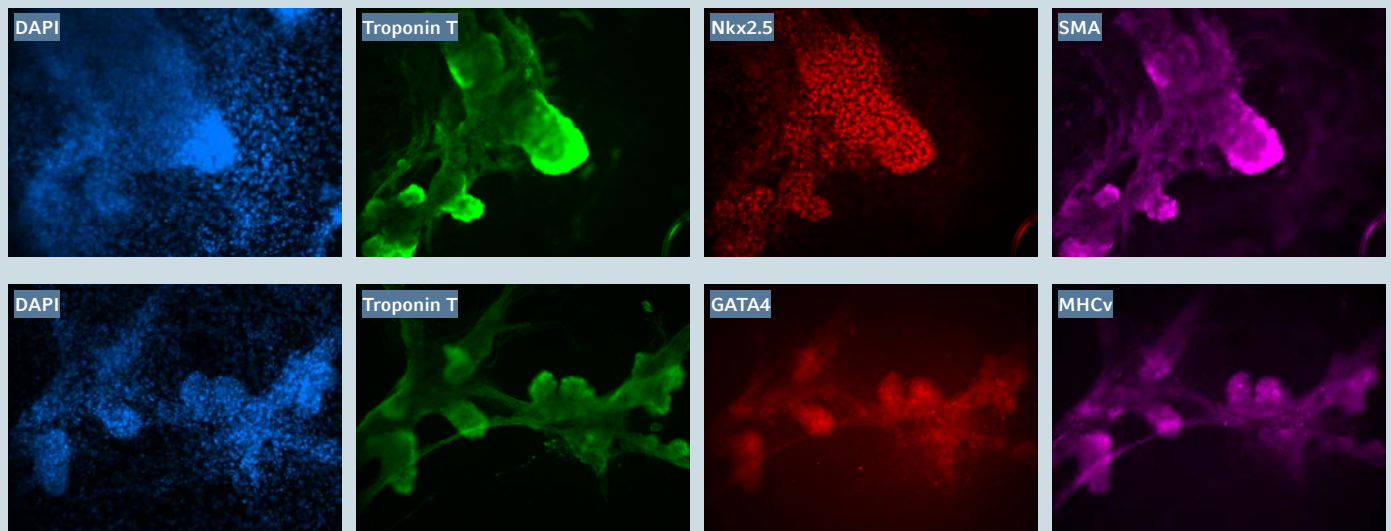
the final differentiation into presumptive motor neurons (Figure 9, lower panel). After 7 days in the final media cocktail the cells organized a fine network of process which were robustly Neurofilament H+, with some remaining TUJ1+ appearance and few GFAP+ glial cells. Many of these cells expressing TUJ1, also expressed canonical markers of motor neurons, such as HB9, and Islet-1 (data not shown). Together these results clearly showed that hypoxic incubation can be used to derive cells of neuroectodermal lineage.



**Figure 9:** Differentiation into presumptive motor neuron precursor cells in low oxygen environment. Immunostaining of PM1Ep5-ERCv2p8 iPSC cells using multi-step motor neuron protocol.

In addition, we were able to first differentiate the cells into definitive endoderm using Wnt pathway activator CHIR99021 using an optimized two step procedure (data not shown), and then finally develop them into cardiomyocytes after Wnt inhibition using wntC-59 (Figure 10). After 10-11 days in culture, while many cells died from the lack of insulin, the surviving cells organized into small patches

and clumps. These “nodes” were highly three dimensional and expressed Troponin T, Nkx2.5 and Smooth muscle actin (SMA). Further, we also observed robust expression of Troponin T along with GATA4 and MHCv. Together this indicates the successful development of the cardiomyocyte lineage in an hypoxic environment.



**Figure 10:** Differentiation into cardiomyocytes in low oxygen environment. Immunostaining of PM1Ep5-ERCv2p9 iPSC cells grown and differentiated using two step cardiomyocyte generation protocol.

## Conclusion

Using the techniques described herein we successfully demonstrated the robust reprogramming of human foreskin fibroblasts in low oxygen tension (5 %) in the CellXpert C170i CO<sub>2</sub> incubator (Figure 11). The “PM1eP5-ERCv2” iPSC line was characterized as karyotypically normal and found to express expected markers of pluripotency Oct4+/SSEA4+ (but not SSEA1-) and Nanog+/Lin28+/Tra-1-60+. Further, the low O<sub>2</sub> environment supported, the cells were able to reliable differentiation of iPSC cells into various stages of ectoderm (not shown), neuronal stem cells, presumptive motor neuron precursors, and finally motor neurons. As well, the hypoxic atmosphere inside the incubator supported the differentiation of the same iPSC line into cardiomyocytes, an endodermal derivative, using standard protocols ([15, 16]).



**Figure 11:** The CellXpert C170i CO<sub>2</sub> Incubator—[Learn more.](#)

## Literature

- [1] Takahashi, K., et al., Induction of pluripotent stem cells from adult human fibroblasts by defined factors. *Cell*, 2007. 131(5): p. 861-72.
- [2] Yu, J., et al., Induced pluripotent stem cell lines derived from human somatic cells. *Science*, 2007. 318(5858): p. 1917-20.
- [3] Carey, B.W., et al., Reprogramming of murine and human somatic cells using a single polycistronic vector. *Proc Natl Acad Sci U S A*, 2009. 106(1): p. 157-62.
- [4] Fontes, A., et al., Generation of human-induced pluripotent stem cells (hiPSCs) using episomal vectors on defined Essential 8 Medium conditions. *Methods Mol Biol*, 2013. 997: p. 57-72.
- [5] Li, D., et al., Optimized Approaches for Generation of Integration-free iPSCs from Human Urine-Derived Cells with Small Molecules and Autologous Feeder. *Stem Cell Reports*, 2016. 6(5): p. 717-728.
- [6] Onder, T.T., et al., Chromatin-modifying enzymes as modulators of reprogramming. *Nature*, 2012. 483(7391): p. 598-602.
- [7] Questa, M., et al., Generation of iPSC line iPSC-FH2.1 in hypoxic conditions from human foreskin fibroblasts. *Stem Cell Res*, 2016. 16(2): p. 300-3.
- [8] Shimada, H., et al., Accelerated generation of human induced pluripotent stem cells with retroviral transduction and chemical inhibitors under physiological hypoxia. *Biochem Biophys Res Commun*, 2012. 417(2): p. 659-64.
- [9] Yoshida, Y., et al., Hypoxia enhances the generation of induced pluripotent stem cells. *Cell Stem Cell*, 2009. 5(3): p. 237-41.
- [10] Yu, J., et al., Efficient feeder-free episomal reprogramming with small molecules. *PLoS One*, 2011. 6(3): p. e17557.
- [11] Yu, J., et al., Human induced pluripotent stem cells free of vector and transgene sequences. *Science*, 2009. 324(5928): p. 797-801.
- [12] Ren, C.P., et al., Establishment of human embryonic stem cell line stably expressing Epstein-Barr virus-encoded nuclear antigen 1. *Acta Biochim Biophys Sin (Shanghai)*, 2005. 37(1): p. 68-73.
- [13] Chen, G., et al., Chemically defined conditions for human iPSC derivation and culture. *Nat Methods*, 2011. 8(5): p. 424-9.
- [14] Kuo, H.H., et al., Negligible-Cost and Weekend-Free Chemically Defined Human iPSC Culture. *Stem Cell Reports*, 2020. 14(2): p. 256-270.
- [15] Bianchi, F., et al., Rapid and efficient differentiation of functional motor neurons from human iPSC for neural injury modelling. *Stem Cell Res*, 2018. 32: p. 126-134.
- [16] Francis, N.L., et al., Peptide-Based Scaffolds for the Culture and Transplantation of Human Dopaminergic Neurons. *Tissue Eng Part A*, 2019.
- [17] Burridge, P.W., A. Holmstrom, and J.C. Wu, Chemically Defined Culture and Cardiomyocyte Differentiation of Human Pluripotent Stem Cells. *Curr Protoc Hum Genet*, 2015. 87: p. 21 3 1-21 3 15.
- [18] Burridge, P.W., et al., Chemically defined generation of human cardiomyocytes. *Nat Methods*, 2014. 11(8): p. 855-60.
- [19] Weber, K., et al., A multicolor panel of novel lentiviral "gene ontology" (LeGO) vectors for functional gene analysis. *Mol Ther*, 2008. 16(4): p. 698-706.
- [20] Meng, X., et al., Efficient reprogramming of human cord blood CD34+ cells into induced pluripotent stem cells with OCT4 and SOX2 alone. *Mol Ther*, 2012. 20(2): p. 408-16.
- [21] Yu, W., et al., Catalytic site remodelling of the DOT1L methyltransferase by selective inhibitors. *Nat Commun*, 2012. 3: p. 1288.
- [22] Bettger, W.J., et al., Rapid clonal growth and serial passage of human diploid fibroblasts in a lipid-enriched synthetic medium supplemented with epidermal growth factor, insulin, and dexamethasone. *Proc Natl Acad Sci U S A*, 1981. 78(9): p. 5588-92.
- [23] Bettger, W.J. and R.G. Ham, Effects of non-steroidal anti-inflammatory agents and antioxidants on the clonal growth of human diploid fibroblasts. *Prog Lipid Res*, 1981. 20: p. 265-8.
- [24] Walthall, B.J. and R.G. Ham, Multiplication of human diploid fibroblasts in a synthetic medium supplemented with EGF, insulin, and dexamethasone. *Exp Cell Res*, 1981. 134(2): p. 303-11.
- [25] Nahm, W.K., L. Zhou, and V. Falanga, Sustained ability for fibroblast outgrowth from stored neonatal foreskin: a model for studying mechanisms of fibroblast outgrowth. *J Dermatol Sci*, 2002. 28(2): p. 152-8.
- [26] Toma, J.G., et al., Isolation and characterization of multipotent skin-derived precursors from human skin. *Stem Cells*, 2005. 23(6): p. 727-37.

- [27] Byrne, J.A., H.N. Nguyen, and R.A. Reijo Pera, Enhanced generation of induced pluripotent stem cells from a subpopulation of human fibroblasts. *PLoS One*, 2009. 4(9): p. e7118.
- [28] Okita, K., et al., A more efficient method to generate integration-free human iPS cells. *Nat Methods*, 2011. 8(5): p. 409-12.
- [29] Li, F., et al., Apoptotic caspases regulate induction of iPSCs from human fibroblasts. *Cell Stem Cell*, 2010. 7(4): p. 508-20.
- [30] Cotten, M., et al., Lipopolysaccharide is a frequent contaminant of plasmid DNA preparations and can be toxic to primary human cells in the presence of adenovirus. *Gene Ther*, 1994. 1(4): p. 239-46.
- [31] Ma, R., et al., Removing endotoxin from plasmid samples by Triton X-114 isothermal extraction. *Anal Biochem*, 2012. 424(2): p. 124-6.
- [32] Appendix 1: Materials and Methods. For supplementary information, contact your local Eppendorf partner.

**Your local distributor: [www.eppendorf.com/contact](http://www.eppendorf.com/contact)**  
Eppendorf AG · Barkhausenweg 1 · 22339 Hamburg · Germany  
[eppendorf@eppendorf.com](mailto:eppendorf@eppendorf.com) · [www.eppendorf.com](http://www.eppendorf.com)

[www.eppendorf.com](http://www.eppendorf.com)