

Feeder-free derivation of induced pluripotent stem cells from multiple somatic cell types in the defined and low-protein TeSR™-E7™ medium

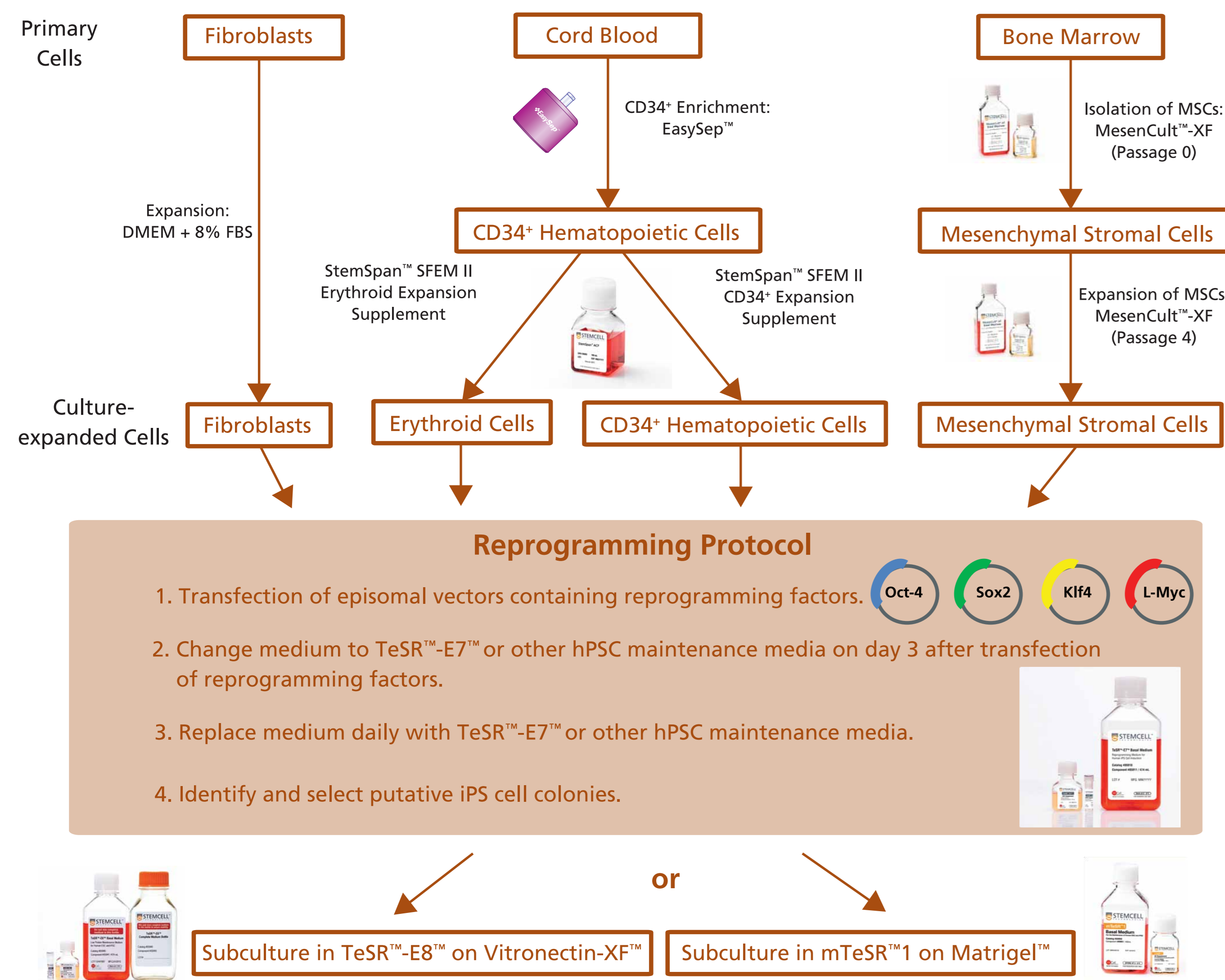
Wing Y. Chang¹, Arwen Hunter¹, Alvin Ng¹, Matthew Wong¹, Susan De Jong¹, Chris Duronio¹, Irene Yu¹, Bert Wognum¹, Ravenska Wagey¹, Carrie Peters¹, Karina McQueen¹, Maureen Fairhurst¹, Erik Hadley¹, Jennifer Antonchuk¹, Terry E. Thomas^{1,2}, Allen C. Eaves^{1,2}, and Sharon A. Louis¹
¹STEMCELL Technologies Inc., Vancouver, Canada ²Terry Fox Laboratory, BC Cancer Agency, Vancouver, B.C., Canada

Introduction

Human induced pluripotent stem (iPS) cells have been derived from a wide range of somatic cell types by the process known as reprogramming. Due to their accessibility, the two most common tissues used for reprogramming are skin (e.g. dermal fibroblasts) and blood (e.g. mononuclear cell fraction, CD34⁺ cells, or specific blood cell types). In addition to identifying a suitable tissue source, there is a general trend towards reprogramming methods that will generate more clinically relevant iPS cells, such as the use of vectors that do not genetically alter the cell, as well as cell culture systems that are free of xenogeneic material and contain defined components. Towards this goal, we recently released TeSR™-E7™ Medium for reprogramming, which is a low-protein, xeno-free, defined medium for reprogramming without the use of feeder cells. This study demonstrates the efficient generation and in depth characterization of iPS cells derived in TeSR™-E7™ Medium from multiple human fibroblast sources, and further extends the use of TeSR™-E7™ Medium for feeder-free reprogramming of other somatic cell types, including bone marrow-derived mesenchymal stromal cells (MSCs), CD34⁺ hematopoietic cells, and culture-expanded erythroid cells. From these cell sources we demonstrate successful establishment of iPS cell lines which can be further cultured in human pluripotent stem cell (hPSC) maintenance media such as mTeSR™1 or TeSR™-E8™. Further confirmation of iPS cell identity in cell lines generated by reprogramming of dermal fibroblasts, includes pluripotency marker expression and in vitro differentiation to cells of the three germ layers. Overall, this study demonstrates the robustness TeSR™-E7™ Medium for reprogramming multiple human somatic cell types, including commonly used tissues such as skin and blood.

Methods

FIGURE 1. Schematic of reprogramming systems tested with TeSR™-E7™



Prior to reprogramming, cells from skin fibroblasts and hematopoietic tissues were isolated and expanded in their specific expansion medium. Adult normal human dermal fibroblasts (NHDF), neonatal fibroblasts (BJ), and fetal fibroblasts (D551) were cultured in fibroblast medium (DMEM + 8% FBS). MSCs were isolated from bone marrow and expanded in MesenCult™-XF Medium. Hematopoietic cells were isolated from cord blood using EasySep™ Human Cord Blood CD34 Positive Selection Kit and expanded in StemSpan™ SFEM II Medium supplemented with either CD34⁺ Expansion Supplement or Erythroid Expansion Supplement. To initiate reprogramming, cells were transfected with episomal vectors containing the reprogramming factors Oct-4, Sox2, Klf4 and L-Myc. Cells were then seeded onto a 6-well plate pre-coated with Matrigel™ and cultured for 3 days in their respective expansion media. After three days, media were replaced with TeSR™-E7™ Medium, which was used for the remainder of the induction phase of reprogramming (day 3 - 28), with daily media changes. Putative iPS cell colonies were typically scored on day 28. iPS cell colonies were manually selected and subcultured either in mTeSR™1 on Matrigel™ or in TeSR™-E8™ Medium on Vitronectin-XF™ for further characterization.

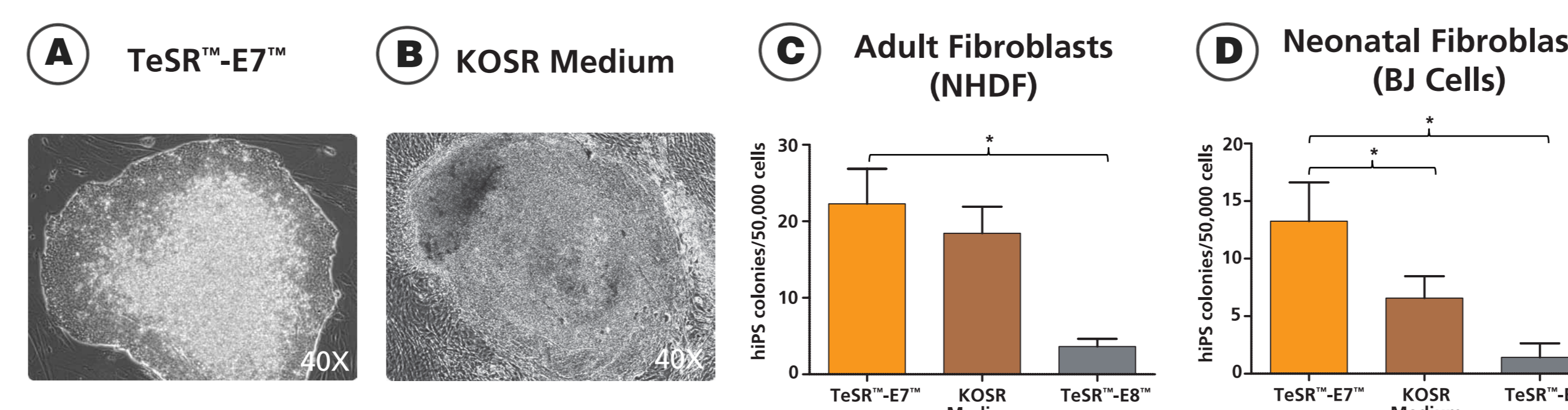
Results

TABLE 1. Fibroblast reprogramming efficiencies in TeSR™-E7™ feeder-free conditions

Somatic Cell Type	# of Experiments	# of Input Cells	Avg. # of iPS Cell Colonies	*Reprogramming Efficiency (%)
Fetal Fibroblasts (D551)	1	5 x 10 ⁴	7	0.015
Neonatal Fibroblasts (BJ)	6	5 x 10 ⁴	13	0.026 ± 0.006
Adult Normal Human Dermal Fibroblasts (NHDF)	6	5 x 10 ⁴	22	0.044 ± 0.009

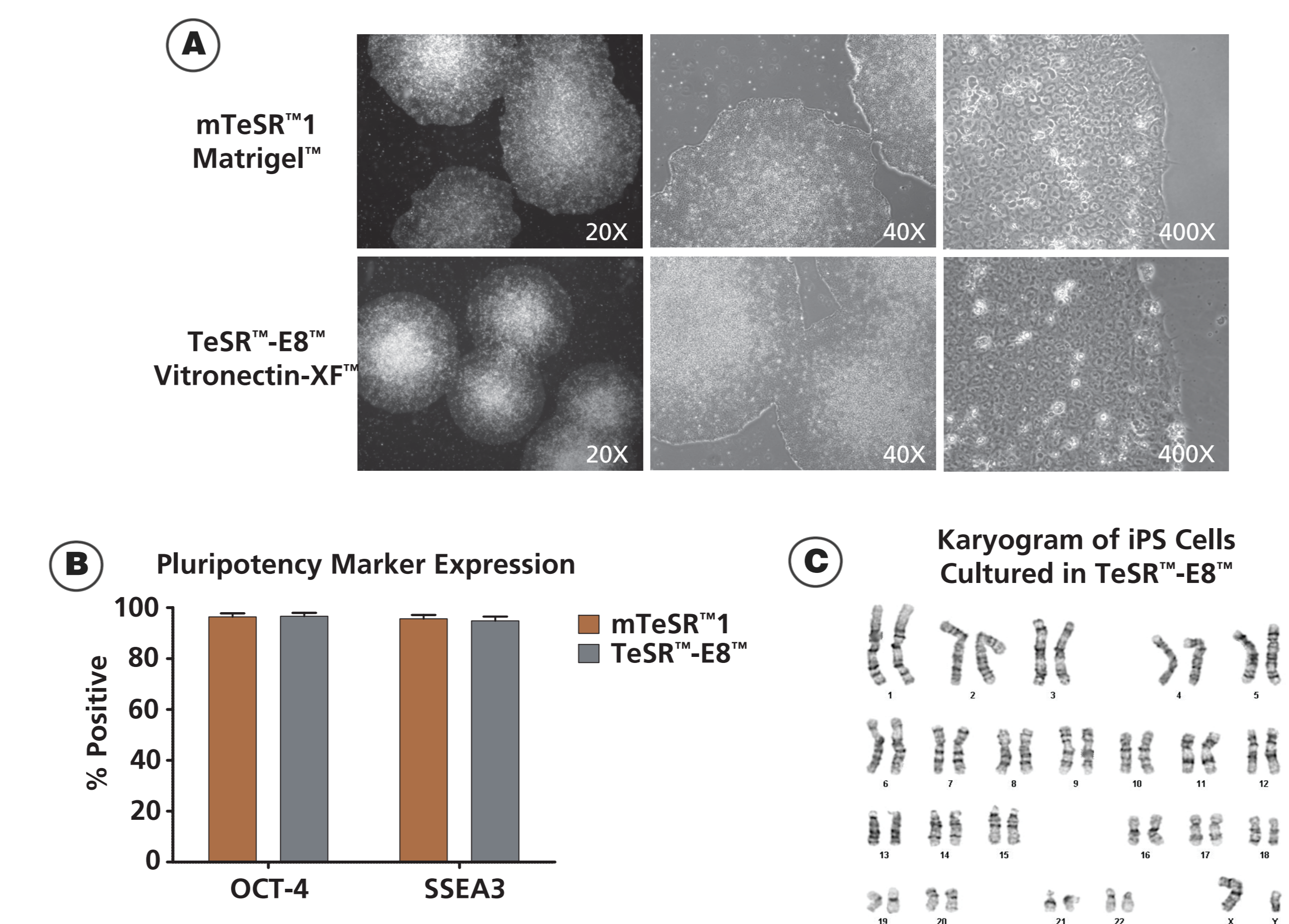
*mean ± SE

FIGURE 2. TeSR™-E7™ yields higher quality iPS cell colonies and improved reprogramming efficiencies compared to standard pluripotent maintenance media



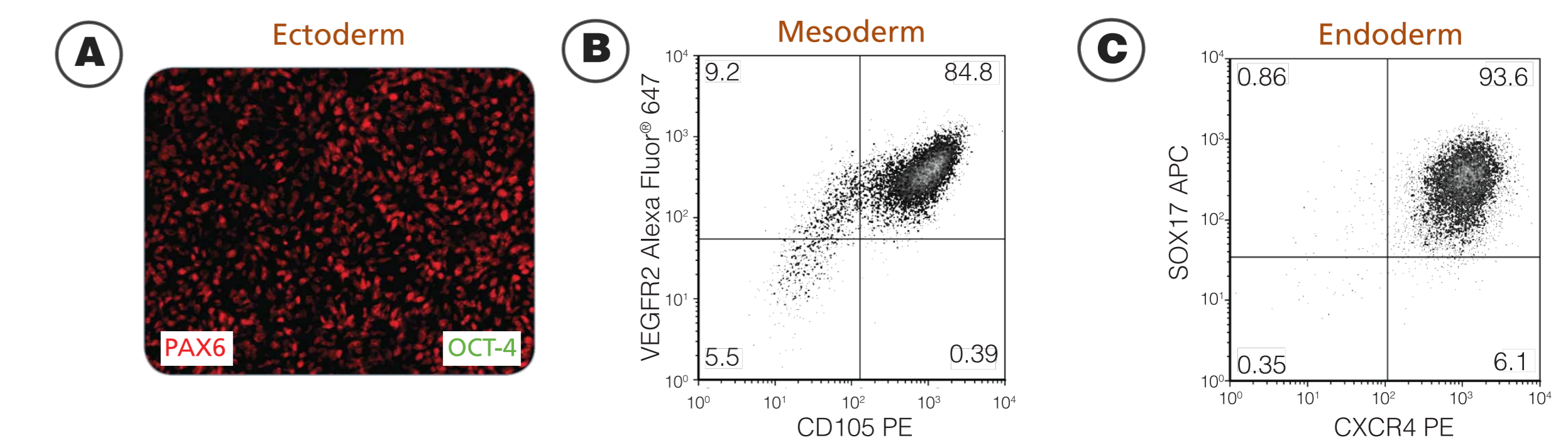
(A - B) Reprogramming of human fibroblasts in TeSR™-E7™ (A) typically yields a greater proportion of high quality iPS cell colonies with defined borders and results in less overgrowth of background fibroblasts compared to KOSR medium (B; DMEM/F12, 20% KnockOut Serum Replacement, bFGF). (C) Reprogramming efficiency for adult fibroblasts in TeSR™-E7™ Medium was comparable to that in KOSR medium, and 6.1-fold higher than in TeSR™-E8™ Medium. (D) For neonatal fibroblasts, reprogramming efficiency in TeSR™-E7™ Medium was 2-fold and 9.4-fold higher than in KOSR medium and TeSR™-E8™ Medium, respectively. Data represents average ± SEM of 6 independent experiments *p<0.05.

FIGURE 3. iPS cells derived in TeSR™-E7™ Medium can be readily subcultured in mTeSR™1 or TeSR™-E8™ Media, expressed pluripotent markers, and exhibited normal karyotype



(A) iPS cell lines derived from dermal fibroblasts in TeSR™-E7™ and subcultured in mTeSR™1 on Matrigel™ or TeSR™-E8™ on Vitronectin-XF™ displayed typical iPS-like morphology with defined borders and high nuclear to cytoplasmic ratio. (B) iPS cell lines derived in TeSR™-E7™ and cultured in either mTeSR™1 or TeSR™-E8™ showed high expression of pluripotency markers SSEA-3 and Oct-4 at passage 2. (C) Karyotype analyses of TeSR™-E7™-derived iPS cells at passage 6 showed no chromosomal anomalies.

FIGURE 4. iPS cells derived in TeSR™-E7™ Medium demonstrated the capacity to differentiate in vitro to cells of the three germ layers



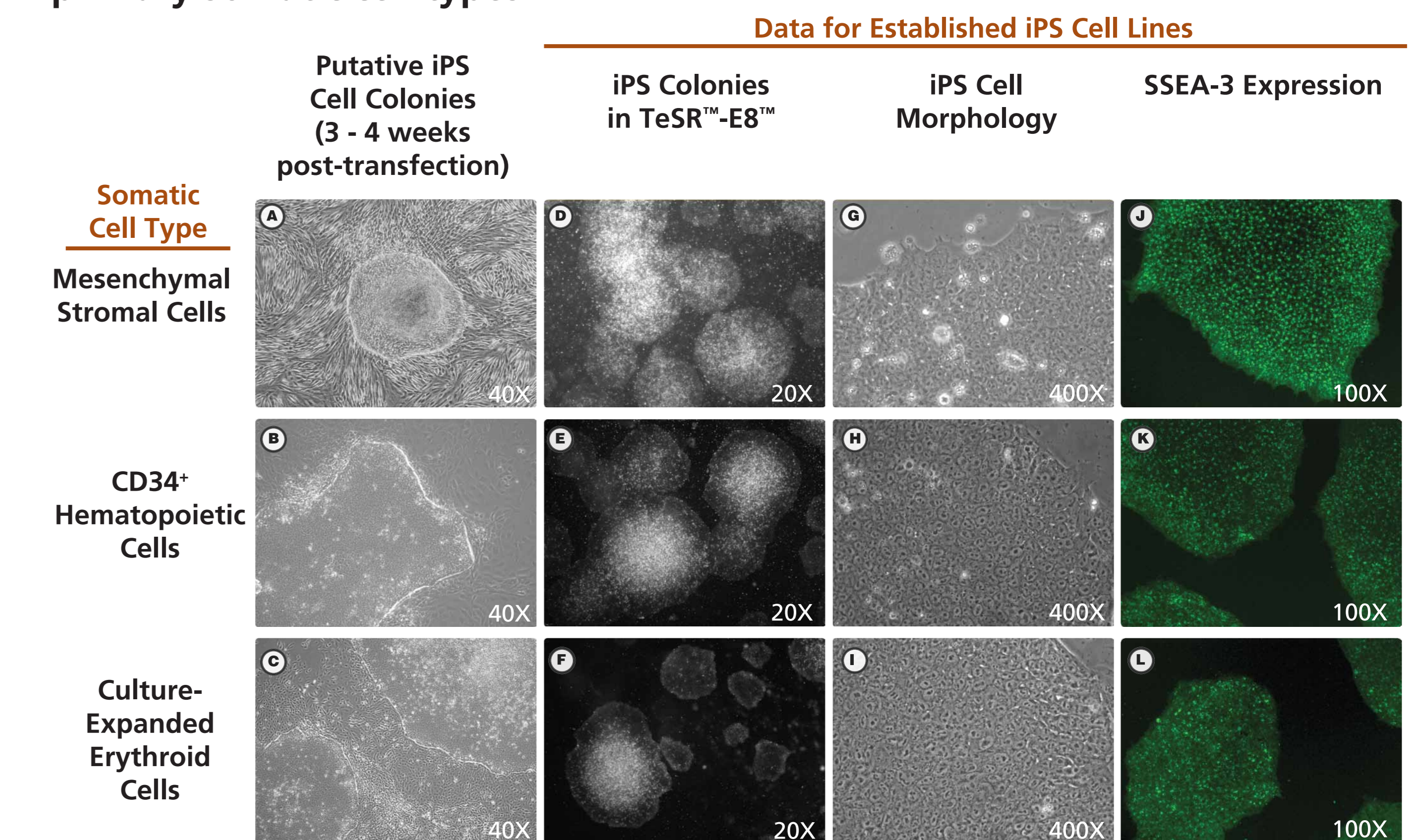
To demonstrate pluripotency, iPS cell lines derived from fibroblasts in TeSR™-E7™ were differentiated to cells of the three germ layers using the STEMdiff™ systems. (A) iPS cells differentiated using STEMdiff™ Neural Induction Medium generated greater than 90% neural progenitor cells expressing PAX6. (B) Mesoderm potential was demonstrated by differentiation of iPS cells to endothelial cells using STEMdiff™ APEL™ Medium supplemented with cytokines. This generated a population of cells with greater than 80% CD105 and VEGFR2 expression. (C) iPS cells differentiated using STEMdiff™ Definitive Endoderm kit exhibited greater than 90% SOX17 and CXCR4-positive cells.

TABLE 2. Reprogramming efficiencies with other somatic cell types in TeSR™-E7™ feeder-free conditions

Somatic Cell Type	# of Input Cells	# of iPS Colonies	*Reprogramming Efficiency (%)
Mesenchymal Stromal Cells	5 x 10 ⁵	5	0.001
CD34 ⁺ Hematopoietic Cells	5 x 10 ⁵	75	0.015
Culture-Expanded Erythroid Cells	5 x 10 ⁵	35	0.007

*n = 1

FIGURE 5. TeSR™-E7™ Medium supports the reprogramming of multiple primary somatic cell types



Mesenchymal stromal cells (top row), CD34⁺ hematopoietic cells (middle row) and culture-expanded erythroid cells (bottom row) were reprogrammed in TeSR™-E7™ Medium with episomal reprogramming vectors as described in Figure 1. (A - C) Three to four weeks post-transfection, putative iPS cell colonies began to emerge and were identified microscopically. (D - L) iPS cells generated from all three somatic cell sources subcultured in either TeSR™-E8™ or mTeSR™1 (data not shown), exhibited iPS-like colony morphology (D - F), high nuclear to cytoplasmic ratio (G - I), and expressed the pluripotency marker SSEA-3 (J - L).

Conclusions

- Improved fibroblast reprogramming efficiency and iPS cell colony morphology was achieved with TeSR™-E7™ Medium compared to common hPSC maintenance media.
- iPS cells can be generated from multiple somatic cell types in TeSR™-E7™ Medium, such as mesenchymal stromal cells, CD34⁺ hematopoietic cells, and culture-expanded erythroid cells.
- TeSR™-E7™ derived iPS cell colonies are easily subcultured in mTeSR™1 or TeSR™-E8™ Media.
- iPS cells derived in TeSR™-E7™ Medium expressed pluripotency markers and had the capacity to differentiate to cells of the three germ layers.