

A FEEDER-INDEPENDENT CULTURE SYSTEM TO CONVERT AND MAINTAIN HUMAN PLURIPOTENT STEM CELLS IN A NAÏVE-LIKE STATE

Arwen L. Hunter¹, Kimberly A. Snyder¹, Terry E. Thomas¹, Allen C. Eaves^{1,2}, Sharon A. Louis¹ and Vivian M. Lee¹
¹STEMCELL Technologies Inc., Vancouver, BC, Canada ²Terry Fox Laboratory, BC Cancer Agency, Vancouver BC, Canada

Introduction

The cells that make up the developing embryo transition through a number of pluripotent states as they become increasingly lineage-restricted. Naïve human pluripotent stem cells (hPSCs) are representative of the inner cell mass of the pre-implantation embryo, whereas primed hPSCs are representative of the inner cell mass of the early post-implantation embryo and are classically dependent on basic fibroblast growth factor (bFGF). The ability to capture these states in vitro requires specialized culture media and protocols.

A simplified and reproducible culture protocol for naïve-like hPSCs in vitro would provide a platform to study the regulatory pathways of naïve and primed pluripotency. To date, conditions that maintain hPSCs in naïve-like states have depended on the use of feeder cells for long-term expansion and results vary depending on the quality of the feeder cells. RSeT™ Feeder-Free (RSeT™-FF) Medium supports the reversion of primed hPSCs to a naïve-like state and supports their long-term feeder-independent maintenance in a bFGF-free system.

Methods

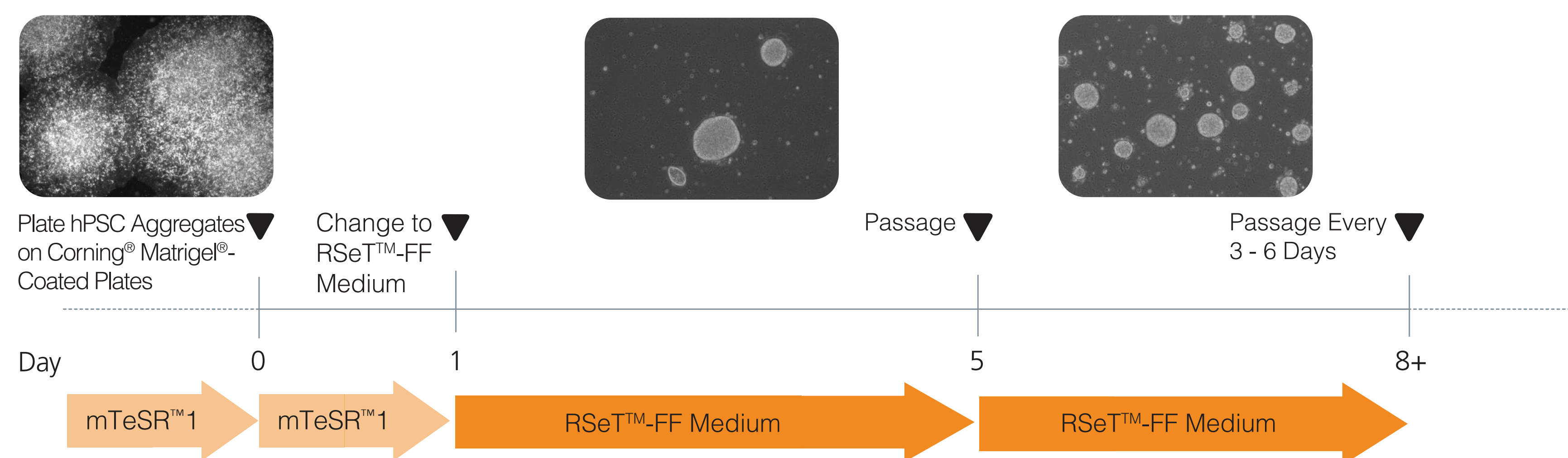


FIGURE 1. RSeT™ Feeder-Free Workflow

Human embryonic stem (ES) and induced pluripotent stem (iPS) cells are reverted to a naïve-like state within one passage in RSeT™-FF Medium. On day 0, seed 500 - 1000 hPSC aggregates per well of a Corning® Matrigel®-coated tissue culture (TC)-treated 12-well plate in mTeSR™1 or TeSR™-E8™. Change medium to RSeT™-FF Medium 24 - 36 hours after seeding. Culture for 4 days in RSeT™-FF Medium and perform full medium changes every other day. Passage cells by incubating cultures with TrypLE™ at 37°C for 8 minutes. Count and plate 20,000 cells/cm² into a non-coated TC-treated plate prepared with RSeT™-FF Medium + 5 µM Y-27632. Perform a full medium change 24 hours after passaging and every other day between passages. Passage cells grown in RSeT™-FF Medium every 3 - 6 days.

Results

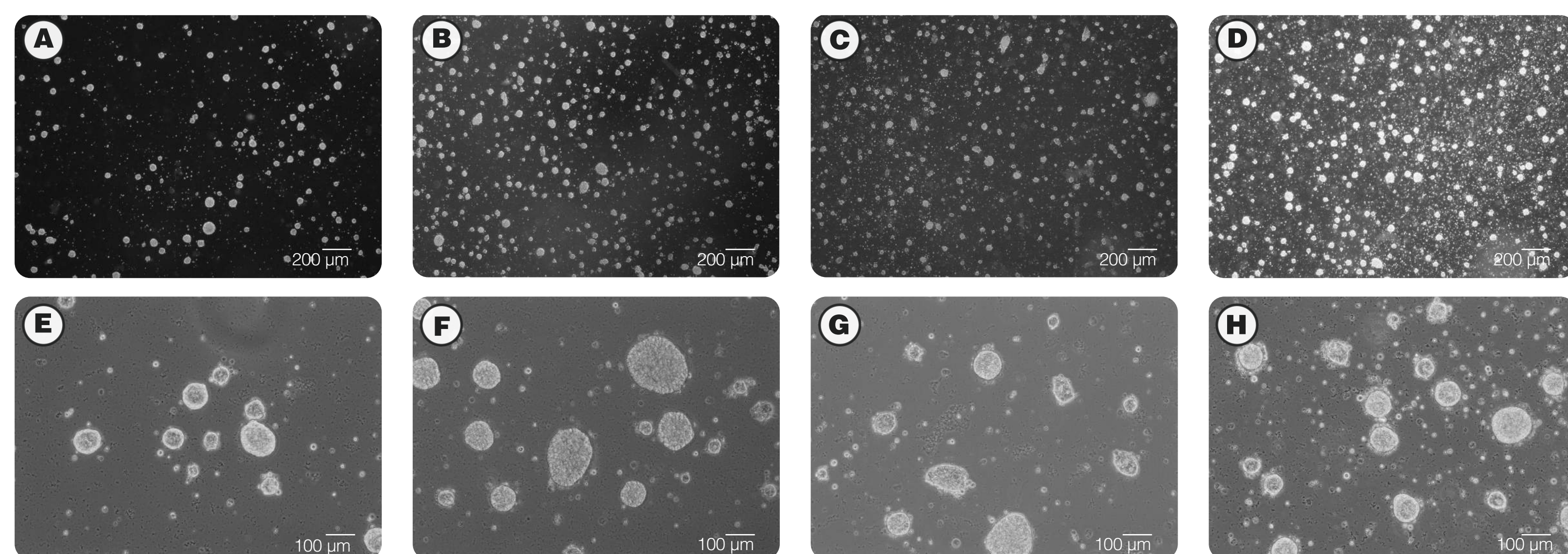


FIGURE 2. hPSCs Cultured in RSeT™ Feeder-Free Medium have a Highly Domed Morphology

Consistent domed morphology of hPSCs cultured in RSeT™-FF Medium in multiple cell lines including: (A, E) H7 hES cells, p10; (B, F) H9 hES cells, p5; (C, G) WLS-1C hiPS cells, p5; and (D, H) STiPS-F016 hiPS cells, p10. These cultures also display very low levels of spontaneous differentiation. Additionally, Corning® Matrigel® supplemented in RSeT™-FF Medium eliminates the need to pre-coat plates.

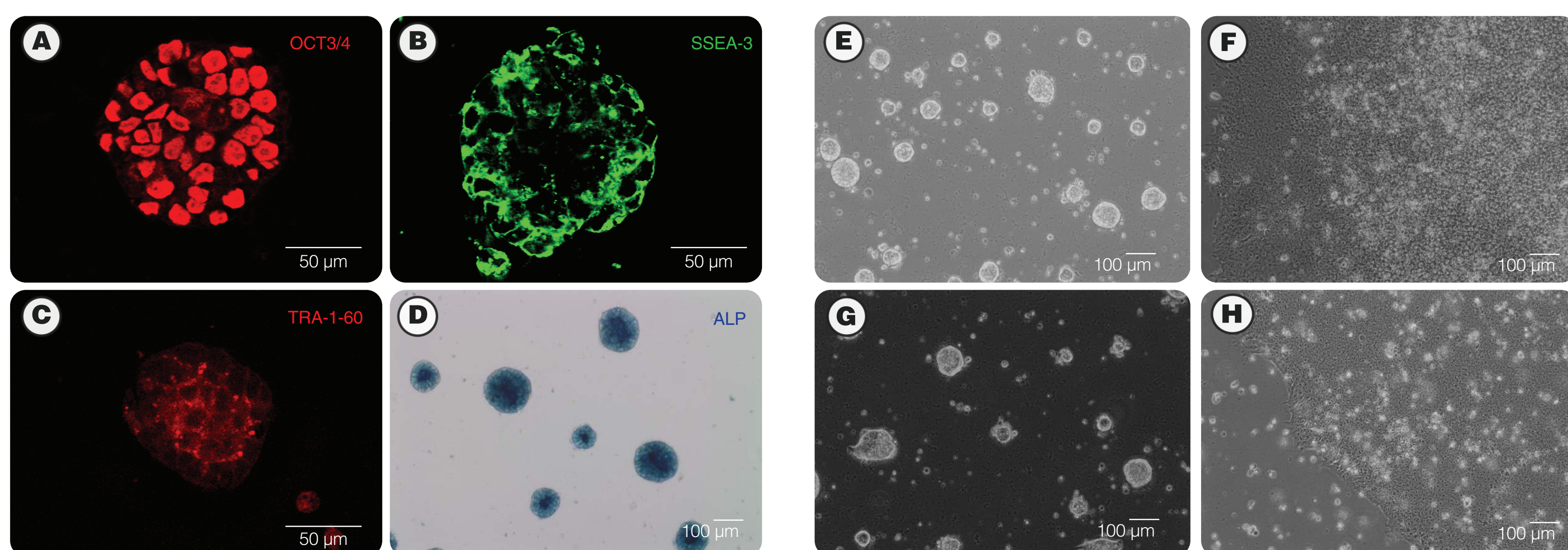


FIGURE 3. hPSCs Cultured in RSeT™-FF Medium Express hPSC Markers and Can Efficiently Transition to a Primed State

Naïve-like hPSCs cultured in RSeT™-FF Medium express (A) OCT3/4 (WLS-1C hiPS cells, p7), (B) SSEA-3 (WLS-1C hiPS cells, p7) (C) TRA-160 (WLS-1C hiPS cells, p7) and (D) stain positive for alkaline phosphatase (ALP) (STiPS-F016 hiPS cell line, p2). Additionally, RSeT™-FF Medium cultures (E: H9 hES cells, G: WLS-1C hiPS cells) can easily transition to a primed state by passaging into either mTeSR™1 (F: H9 hES cells, p1 in mTeSR™1) or TeSR™-E8™ (H: WLS-1C hiPS cells, p2 in TeSR™-E8™).

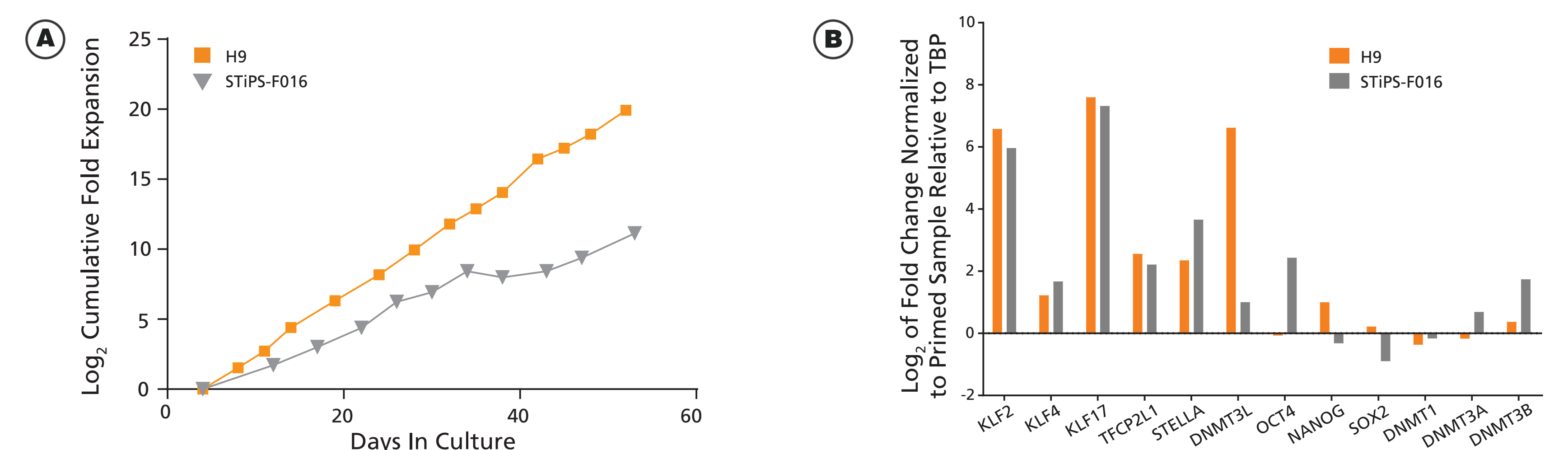


FIGURE 4. hPSCs Cultured in RSeT™-FF Medium Express Genes Associated with the Naïve-Like State

hPSCs maintained in RSeT™-FF Medium (A) display continued expansion and (B) express higher levels of KLF2, KLF4, KLF17, TFPC2L1, STELLA, and DNMT3L compared to primed hPSC controls. Genes associated with pluripotency (OCT4, NANOG, and SOX2) are expressed by naïve-like hPSCs at similar levels compared to primed hPSC controls.



FIGURE 5. hPSCs Maintained in RSeT™-FF Medium Retain a Normal Karyotype After 10 Passages

Representative karyogram results for a (A) H9 hES cells and (B) STiPS-F016 hiPS cells cultured in RSeT™-FF Medium for 10 passages.

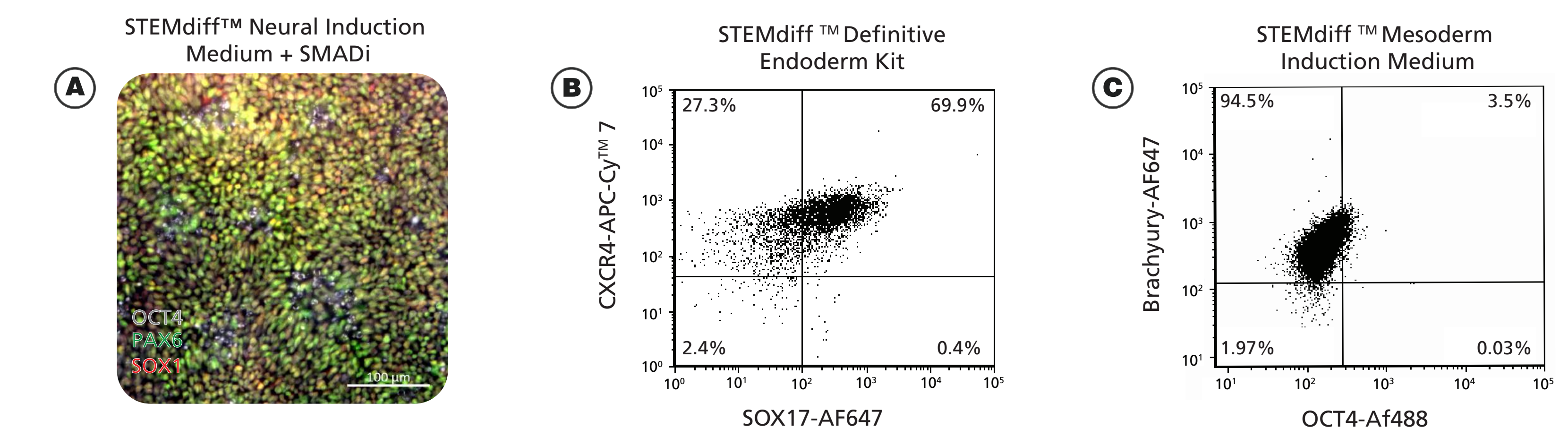


FIGURE 6. hPSCs Maintained in RSeT™-FF Medium are Capable of Direct Trilineage Differentiation Without Re-Priming

(A) Representative image of WLS-1C hiPS cells previously maintained in RSeT™-FF Medium for 11 passages and subjected to neural induction using STEMdiff™ Neural Induction Medium + SMADi monolayer culture. Cultures were processed for immunocytochemistry on day 6 where the majority of cells were positive for neural progenitor markers PAX6 (green) and SOX1 (red), and negative for OCT4 (grey). (B) Representative flow cytometry plot for H9 human ES cells previously maintained in RSeT™-FF Medium and subjected to differentiation using the STEMdiff™ Definitive Endoderm Kit. Cultures were processed for flow cytometry on day 5 where 69.9% of cells were SOX17⁺/CXCR4⁺, indicating differentiation to endodermal lineage. (C) Representative flow cytometry plot for STiPS-F019 hiPS cells previously maintained in RSeT™-FF Medium for 23 passages and subjected to differentiation using STEMdiff™ Mesoderm Induction Medium. Cultures were processed for flow cytometry on day 5 where 94.5% of cells were T⁺/OCT4⁻, indicating efficient differentiation to mesoderm lineage.

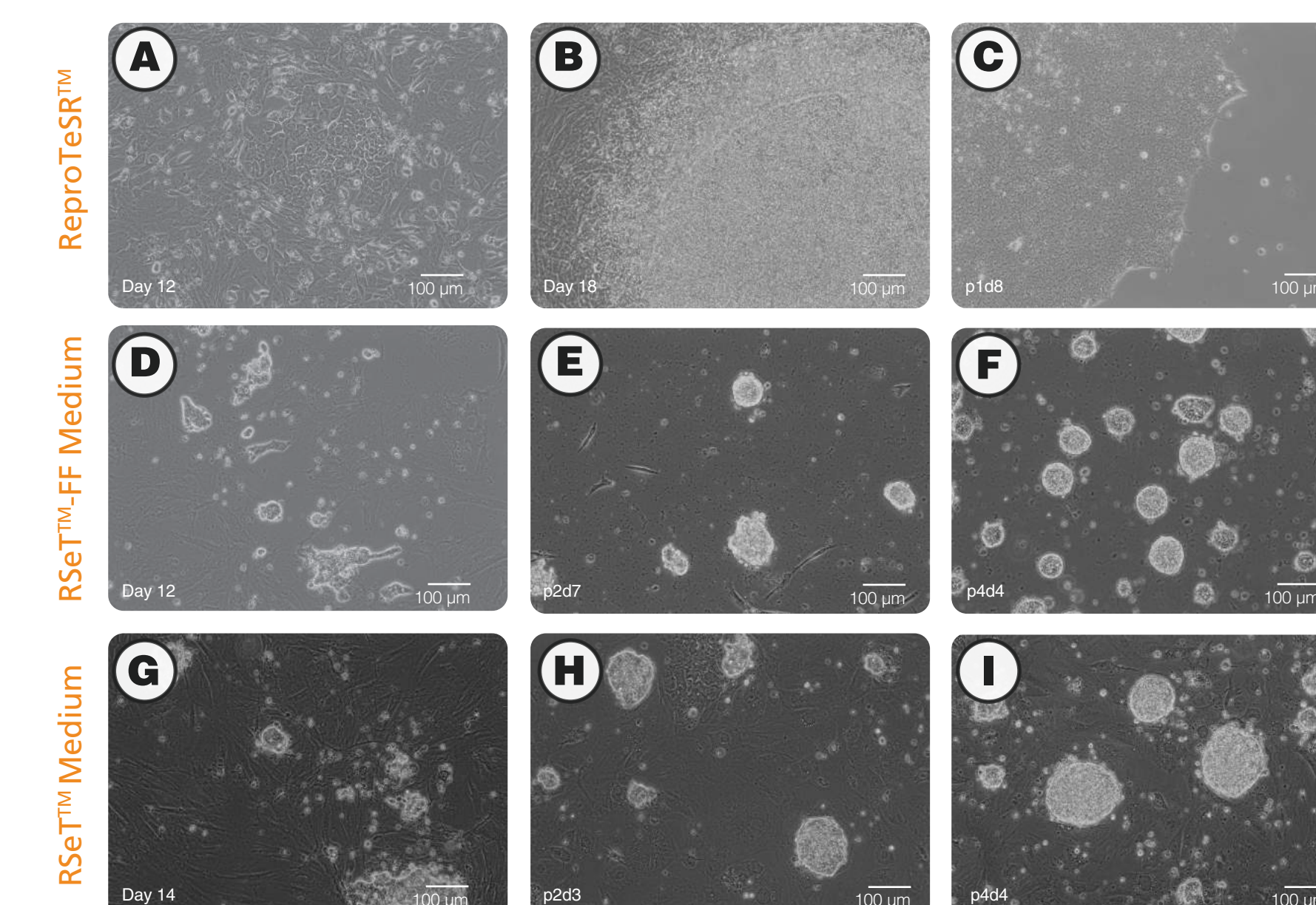


FIGURE 7. Normal Human Dermal Fibroblasts (NHDFs) can be Efficiently Reprogrammed Directly to the Naïve-Like State in Both RSeT™-FF Medium and RSeT™ Medium

Colonies that arise from NHDFs reprogrammed using ReptoRNA™-OKSGM + B18R and either ReptoTeSR™ (A, B), RSeT™-FF Medium (D), or RSeT™ Medium (G). Cultures were passaged on day 12 (RSeT™-FF Medium) or day 14 (RSeT™ Medium) and could be propagated long-term with domed morphology and low levels of differentiation (E, F and H, I; respectively). On day 18, the ReptoTeSR™ control (B) contained colonies large enough to be isolated and propagated in mTeSR™1 (C).

Summary

- RSeT™-FF Medium efficiently reverts primed hPSCs to a naïve-like state with stable domed morphology, naïve gene expression profiles, and low levels of spontaneous differentiation as early as passage 1
- RSeT™-FF Medium maintains hPSCs in a naïve-like state without bFGF or feeder cells in long-term cultures (>p 25)
- Somatic fibroblasts can be reprogrammed directly to the naïve-like state in RSeT™ Medium and RSeT™-FF Medium