

A Simple, Reproducible Method to Generate Red Blood Cells From Human Pluripotent Stem Cells

Crystal Chau¹, Marta A. Walasek¹, Matthew Richardson¹, Christian Barborini¹, Theresa Tai¹, Chelsea Williams¹, Stephen J. Szilvassy¹, Sharon A. Louis¹, Allen C. Eaves^{1,2}, and Albertus W. Wognum¹

¹STEMCELL Technologies Inc., Vancouver, Canada; ²Terry Fox Laboratory, BC Cancer Agency, Vancouver, Canada

INTRODUCTION

Erythroid cells generated from human pluripotent stem cells (hPSCs) offer a potentially unlimited and safe supply of red blood cells (RBCs) for use in disease modeling, drug development, and transfusion medicine. However, obtaining sufficient quantities of RBCs is hampered by technical challenges and the lack of reproducible and robust protocols to differentiate hPSCs into erythroid cells. We have developed a simple two-step, serum- and feeder cell-free culture system (STEMdiff™ Erythroid Kit) to generate erythroid cells from hPSCs. Using only three supplements that combine cytokines and other factors to support optimal cell expansion and differentiation, this method generates a large number of erythroid cells from multiple hPSC lines within 31 days. hPSCs are first differentiated into hematopoietic progenitor cells (HPCs) by specification to mesoderm and hemoendothelial lineage in a monolayer culture system. Next, hPSC-derived HPCs are expanded and differentiated into CD71^{high} GlyA⁺ erythroblasts using erythroid lineage-specific culture conditions. Further culture of these cells results in a > 90% pure population of GlyA⁺ erythroid cells and no cell loss. The cells generated in this system exhibit orthochromatic normoblast morphology and decreased CD71 expression, consistent with erythroid maturation. In summary, we have developed a simple, standardized, and reproducible two-step, serum- and feeder cell-free method for generating large numbers of erythroid cells from multiple hPSC lines for basic and translational research.

METHODS

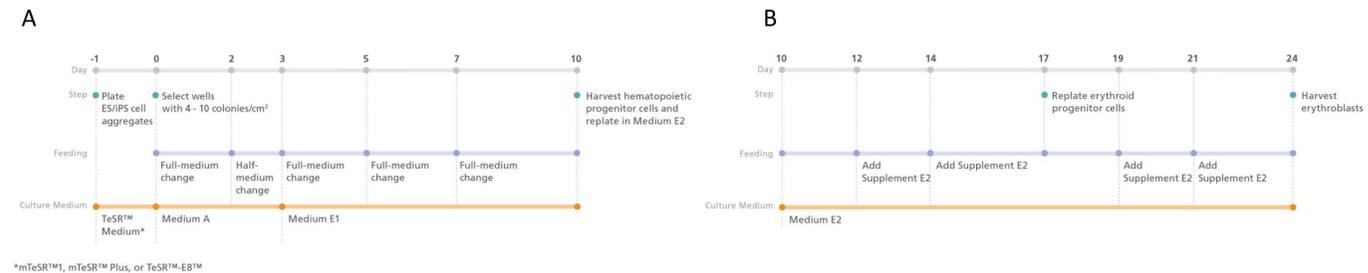


FIGURE 1. Culture Protocol for Generating Erythroblasts from hPSCs

The protocol for the STEMdiff™ Erythroid Kit involves two steps: **(A)** Hematopoietic specification of hPSCs, and **(B)** differentiation of hPSC-derived HPCs into erythroid cells. **(A)** Human embryonic stem (ES) or induced pluripotent stem (iPS) cells were plated as small aggregates (100 - 200 μm diameter) at a density of 15 - 20 aggregates/cm² in mTeSR™1, mTeSR™ Plus, or TeSR™-E8™ medium on Corning® Matrigel®-coated plates, and allowed to attach overnight. The next day, mesoderm differentiation was initiated by changing the medium to Medium A (STEMdiff™ Hematopoietic Basal Medium + STEMdiff™ Hematopoietic Supplement A). After 3 days, the medium was changed to Medium E1 (STEMdiff™ Hematopoietic Basal Medium + STEMdiff™ Erythroid Supplement E1) to induce hemoendothelial and hematopoietic specification. On day 10, HPCs were harvested from the suspension culture and counted. **(B)** Hematopoietic progenitors were replated in Medium E2 (StemSpan™ SFEM II + STEMdiff™ Erythroid Supplement E2) at a density of 4 x 10⁴ cells/mL and the cells were cultured for 14 days to generate erythroblasts. Cells were harvested from the cultures at multiple timepoints, counted, and analyzed by flow cytometry for expression of hematopoietic (CD43 and CD34) and erythroid (CD71 and GlyA) markers. Dead cells were excluded by light-scatter profile and propidium iodide staining. The number of erythroid cells was calculated as the fraction of cells expressing GlyA on their surface. Erythroid identity of the generated cells was additionally assessed by cell morphology and hemoglobin expression. Six different hPSC lines were used in this study: the embryonic stem cell lines H1, H7, and H9, and the induced pluripotent stem cell lines 1C, F016, and B004. All hPSC lines were maintained in feeder-free and serum-free culture conditions prior to differentiation.

RESULTS

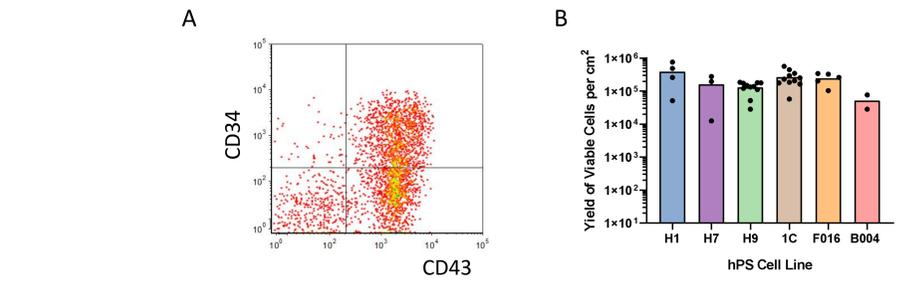


FIGURE 2. hPSCs Differentiate into Hematopoietic Progenitor Cells with High Yields

The hematopoietic specification protocol described in Figure 1A promotes the generation of large numbers of hematopoietic progenitor cells. On day 10, the supernatant cells were harvested and the phenotype and cell yields were calculated **(A&B)**. **(A)** The average frequency of cells expressing the embryonic pan-hematopoietic marker CD43 is 88% (range: 65 - 99%) and the frequency of CD34⁺ cells ranges between 24 - 88%. A representative flow cytometry plot of day 10 hematopoietic progenitors derived from F016 iPS cells is shown. **(B)** The yield of total viable cells per cm² from 6 hPSC lines. The overall average across all cell lines was 2.2 x 10⁵ cells/cm². Black dots depict the results of individual experiments, and each dot is an average of two technical replicates; bars show the mean.

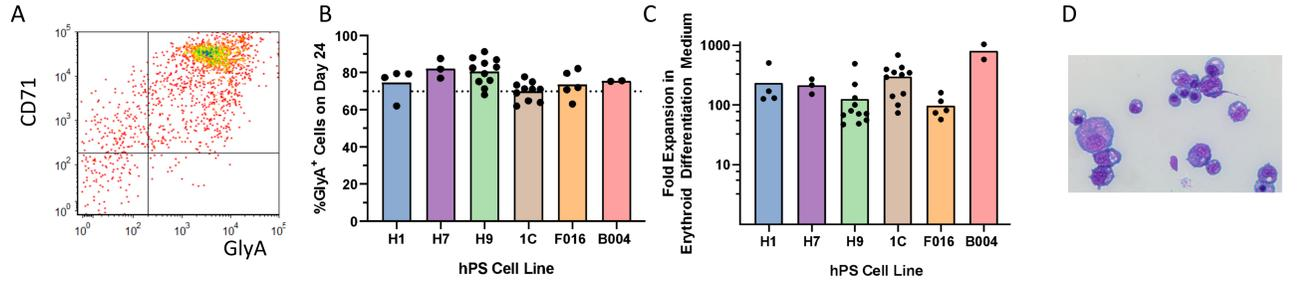


FIGURE 3. hPSC-Derived Hematopoietic Progenitor Cells are Effectively Expanded and Differentiated into Erythroid Cells

hPSC-derived HPCs were differentiated into the erythroid lineage as described in Figure 1B. After 14 days in erythroid culture conditions, the cells were harvested, counted, and analyzed for expression of the erythroid markers CD71 and GlyA **(A-C)**. **(A)** A representative flow cytometry plot of F016 iPS-derived erythroid cells after a total of 24 days of culture. **(B)** The average frequencies of GlyA⁺ erythroid cells on day 24 across 6 hPSC lines. The overall average %GlyA⁺ cells across all cell lines was 76%. Black dots depict the results of individual experiments, and each dot is an average of two technical replicates; bars show the mean. **(C)** Erythroid differentiation conditions supported high expansion of hPSC-derived HPCs with overall average 230-fold expansion across all hPSC lines. Black dots depict the results of individual experiments, and each dot is an average of two technical replicates; bars show the mean. **(D)** Representative image taken at 40X magnification of May-Grunwald-Giemsa stained cells cultured for a total of 24 days (F016 iPS cell line).

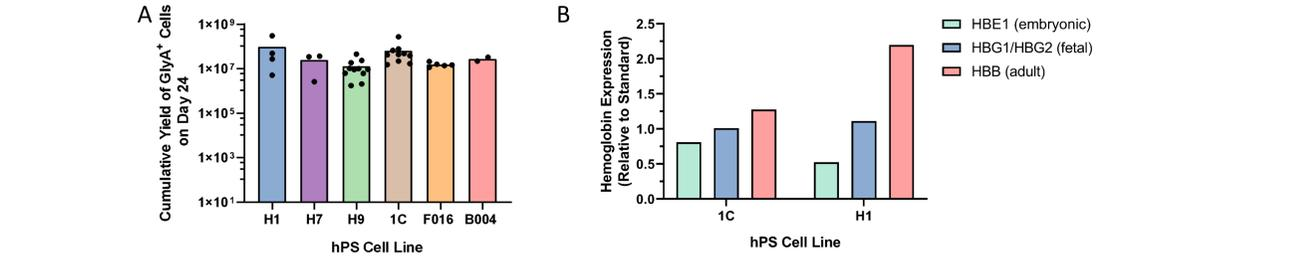


FIGURE 4. Large Numbers of Erythroid Cells Expressing a Mix of Primitive and Definitive Hemoglobin are Generated From hPSC-Derived Hematopoietic Progenitor Cells

(A) Optimal cell differentiation and expansion during both the hematopoietic and erythroid specification steps of the protocol results in the generation of large numbers of erythroid cells after a total of 24 days of culture. The overall average number of generated erythroid cells was 4 x 10⁷ GlyA⁺ cells/cm² across all hPSC lines. Shown are the cumulative yields of GlyA⁺ cells per cm² of the initial culture. Black dots depict the results of individual experiments, and each dot is an average of two technical replicates; bars show the mean. **(B)** Erythroid cells generated on day 24 of the differentiation protocol expressed a mix of 'primitive' (HBE1) and 'definitive' (HBG1/HBG2 and HBB) hemoglobin types, but with adult and fetal hemoglobin expressed at higher levels than embryonic hemoglobin. Representative RT-qPCR data are shown for two cell lines; 1C (iPS) and H1 (ES). Hemoglobin expression was normalized to the expression of housekeeping genes 18S and TBP.

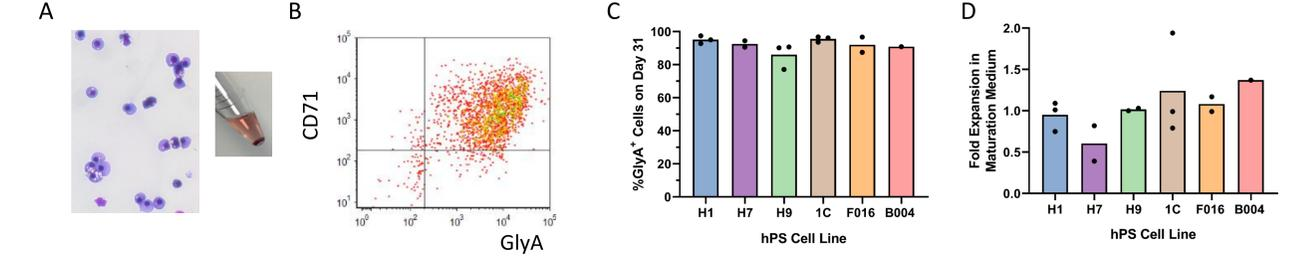


FIGURE 5. hPSC-Derived Erythroblasts Can Mature in Culture

Erythroblasts generated from hPSC-derived HPCs were further cultured for 7 days under maturation conditions with 3 U/mL erythropoietin (EPO) + 3% human serum. After a total of 31 days of culture, the cells were harvested, counted, and assessed for expression of erythroid markers CD71 and GlyA, and changes in cell morphology. **(A)** Differentiated cells were characterized by orthochromatic normoblast morphology (left, 40X magnification), the red color of the cell pellet (right), and **(B)** decreased CD71 expression, which are all consistent with erythroid maturation. **(B&C)** Maturation cultures resulted in a ≥ 90% pure population of GlyA⁺ erythroid cells. **(B)** Representative flow cytometry plot of cells derived from F016 iPS cells. **(C)** Bar graphs summarize the average frequencies of GlyA⁺ erythroid cells on day 31 across all 6 hPSC lines. Black dots depict the results of individual experiments, and each dot is an average of two technical replicates; bars show the mean. **(D)** Notably, no cell loss was observed during the maturation culture. Black dots depict the results of individual experiments, and each dot is an average of two technical replicates; bars show the mean. The overall average cumulative yields of GlyA⁺ cells after 31 days of culture were 9.3 x 10⁷ cells per initial cm² across all cell lines (data not shown).

Summary

- hPSCs can be differentiated into erythroid cells with the STEMdiff™ Erythroid Kit, a simple two-step, feeder cell-free and serum-free culture system. The STEMdiff™ Erythroid Kit will be launched in early 2021.
- Each culture step supports optimal cell expansion and differentiation efficiency and results in high purity and yields of GlyA⁺ erythroid cells from multiple hPSC lines.
- The generated erythroid cells expressed a mix of adult, fetal, and embryonic hemoglobin types.
- Following maturation culture, the cells display orthochromatic normoblast morphology.
- Future work will aim to further increase erythroid cell maturation in vitro.