

# Enabling Efficient Ex Vivo Expansion of Long-term Human Hematopoietic Stem and Progenitor Cells For Cell Therapy

Janet J. Sei, Abigail H. Becker, Navjot Kaur, Mohan C. Vemuri, and David T. Kuninger. Thermo Fisher Scientific, Cell Biology, Frederick MD, USA, 21704.

## ABSTRACT

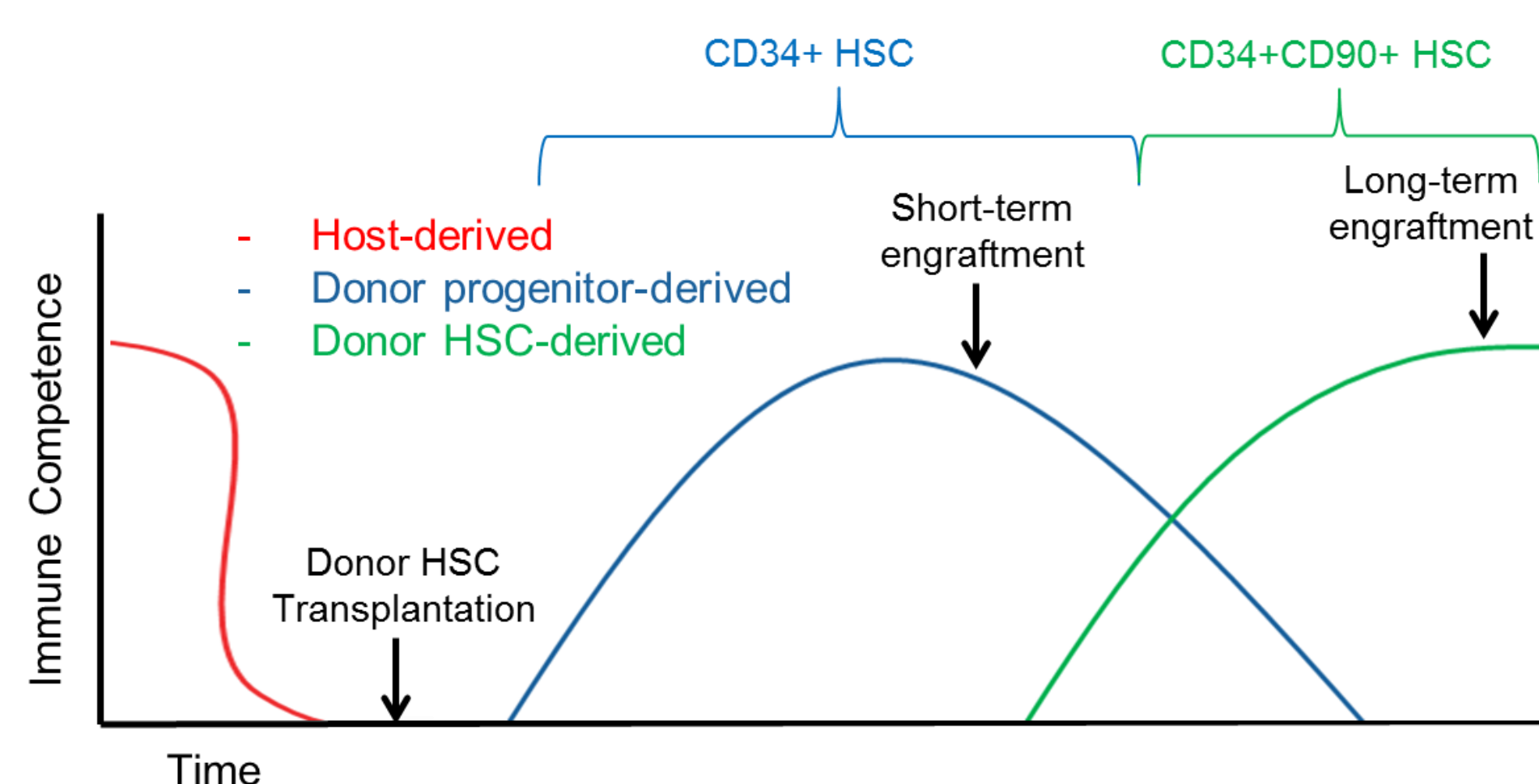
Umbilical-cord blood (UCB) is an important source of hematopoietic stem and progenitor cells (HSPC) for transplantation into patients lacking a suitable HLA-matched donor. However, due to the limited cell dose in each cord blood unit, individuals >60 kg are restricted from the use of UCB-based therapy. Ex vivo expansion of UCB CD34<sup>+</sup> cells is one strategy employed to increase the hematopoietic cell dosage. A major limitation of current systems used for the expansion of HSPC is that ex vivo culture leads to expansion and differentiation at the expense of the most primitive pluripotent long-term stem cells. This has limited the clinical application of ex vivo expanded HSPC, since short-term progenitor cells only provide transient protection, ultimately reducing the long-term positive health outcomes, increasing the duration of hospitalizations, and health care costs per patient. Development of a culture system that expands both short-term and long-term HSPC would facilitate immune protection during the early phase of recovery, and provides a suitable solution for transfusion-independent hematopoiesis. Therefore, we sought to develop a HSPC culture medium that enables the expansion of both long-term and short-term HSPC, while maintaining their functional properties. To this end, we conducted several iterative rounds of Design of Experiments (DOE) involving multifactorial analysis, and mathematical modeling methods. Definitive Screening DOEs allowed us to identify optimal combinations and concentrations of essential media components, small molecules, and growth factors. The performance of candidate HSPC expansion media were evaluated after 7 days of culture, with the following attributes assessed: (1) viability of cells; (2) numbers of total nucleated cells; (3) percentages and numbers of CD34<sup>+</sup> cells; (4) percentages and numbers of CD34<sup>+</sup>CD90<sup>+</sup>CD45RA<sup>-</sup> cells; (5) expression of aldehyde dehydrogenase by expanded CD34<sup>+</sup> cells; and (6) colony-forming unit (CFU) assays. As the transplantation of HSPC in immuno-deficient mice is the gold standard in determining whether the expanded cells are engraftable, we plan to conduct these studies with the lead candidate HSPC expansion medium. Taken together, we seek to highlight our design philosophy in HSPC culture media development. We believe that our efforts are critical for the successful utilization of hematopoietic stem cell transplants in translational cell therapies.

## INTRODUCTION

The goal of this research was to develop a hematopoietic stem and progenitor cell (HSPC) expansion media that fulfills the needs for customers interested in therapies that treat cancer, blood disorders, immunological disorders, in addition to generation of iPSC and cardiomyocyte.



Current systems used for the ex vivo expansion of HSPC result in the expansion and differentiation of CD34<sup>+</sup> cells, at the expense of the most primitive pluripotent long-term stem cells CD34<sup>+</sup>CD90<sup>+</sup>CD45RA<sup>-</sup> cells that provide life-long immunity.



Adapted from Bhattacharya et al.

## MATERIALS AND METHODS

Figure 1. Process Workflow for Expansion of CD34<sup>+</sup> cells

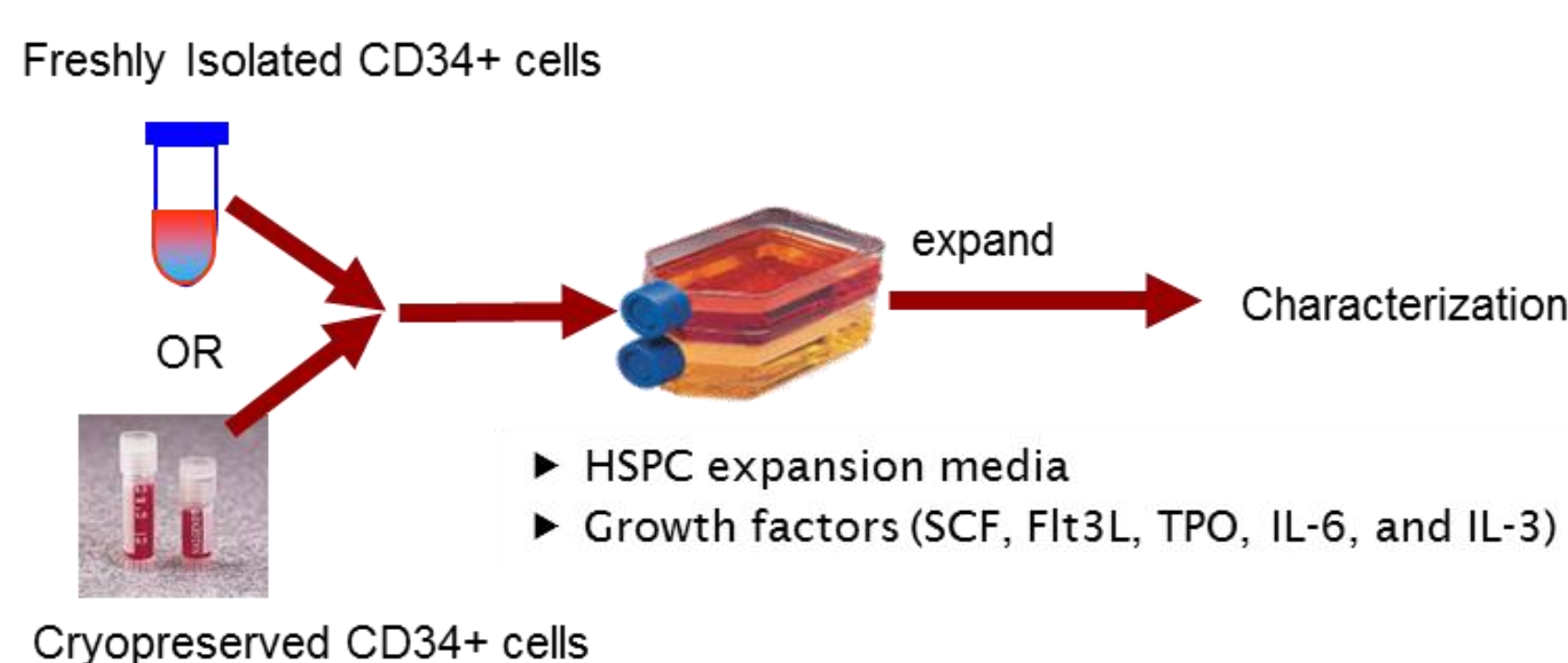


Figure 1: CD34<sup>+</sup> cells were expanded with HSPC expansion media and characterized to determine the TNC, % Viability, phenotype and function.

Figure 2. Experimental Workflow for Development of HSPC Expansion Medium

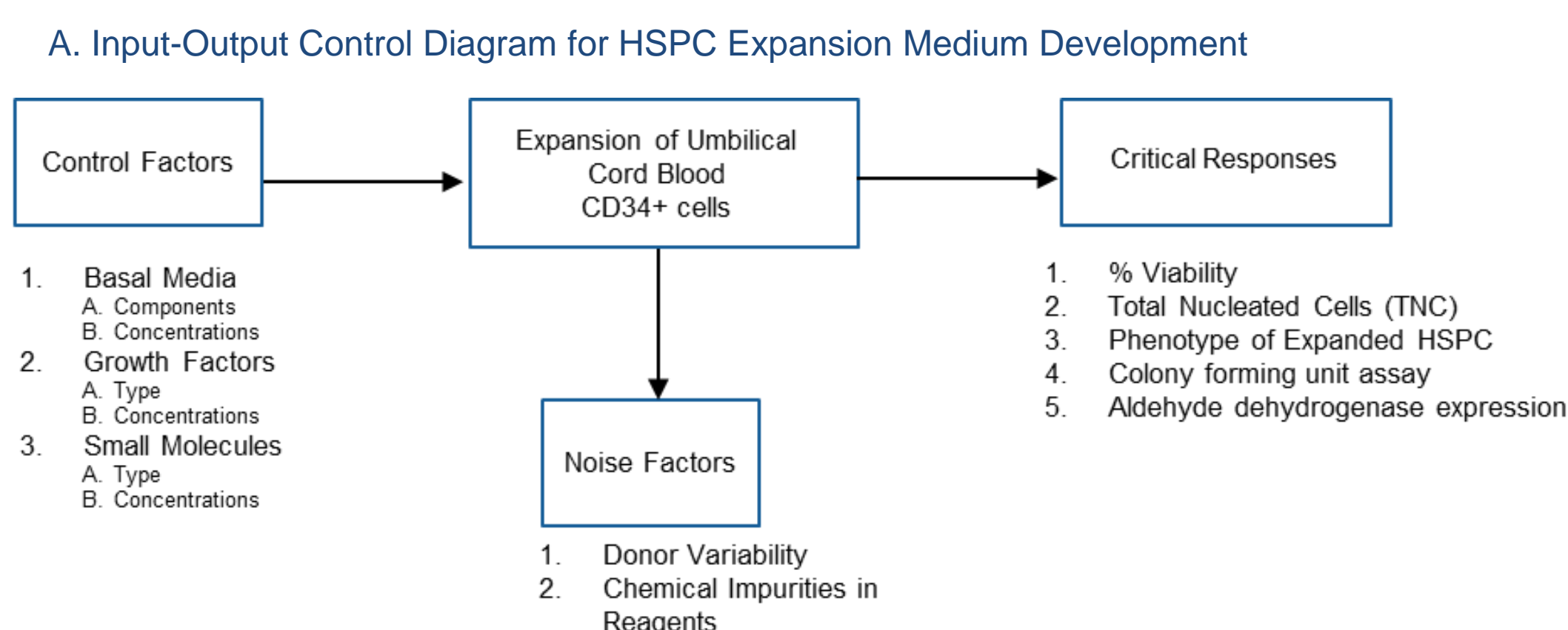


Figure 2: Process Map that outlines input variables that may influence outputs. Illustrates potential Noise Factors that increase variability during medium development.

## RESULTS

Figure 3. Analysis of Existing Solutions in Expansion of Umbilical Cord Blood CD34<sup>+</sup> cells.

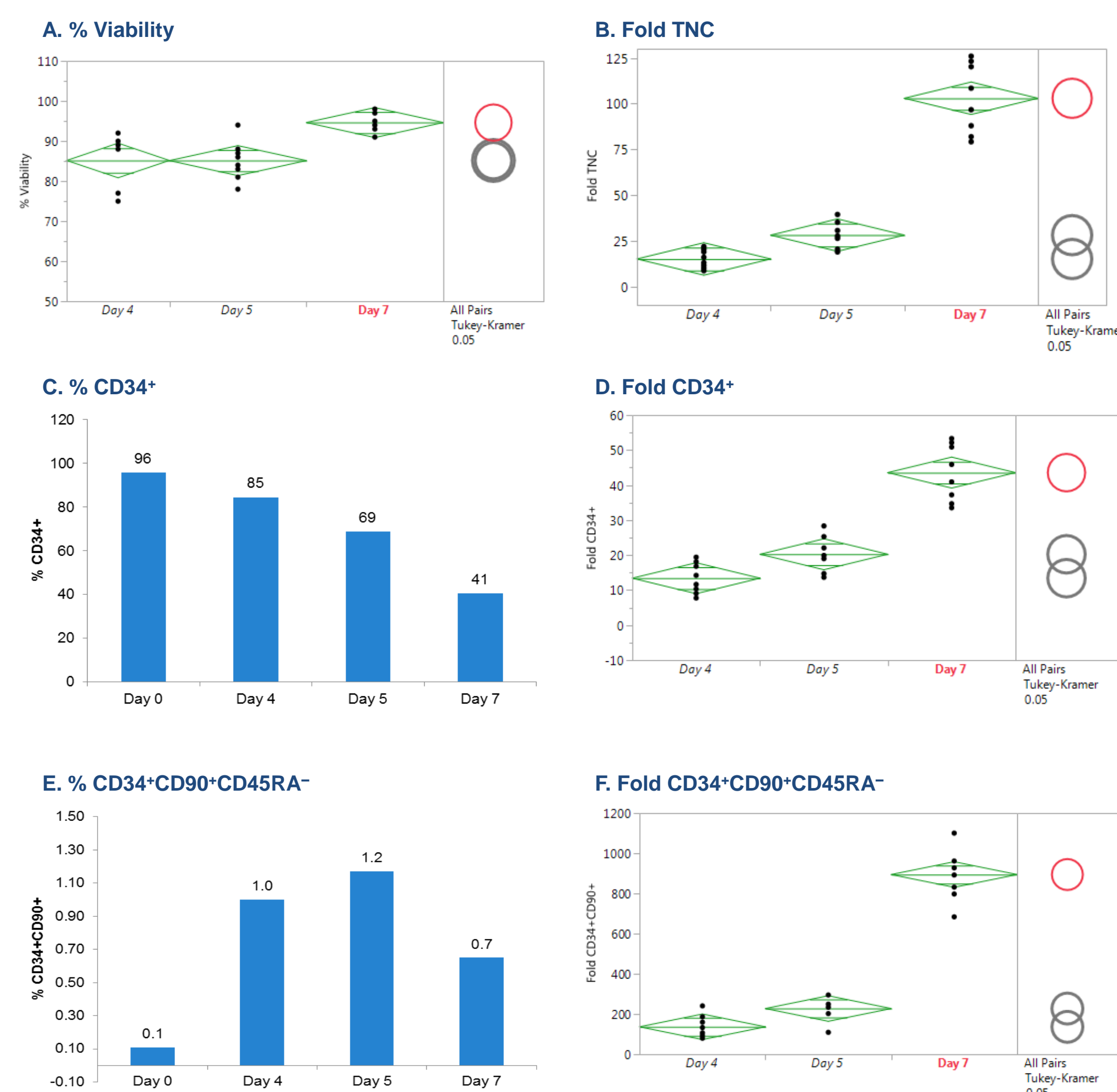


Figure 3: Representative results of current solutions demonstrated that the highest increase in HSPC was observable at 7 days post expansion. Viability of cells is highest at Day 7 (A). While the number of TNC increases with time (B), the % of CD34<sup>+</sup> decreases due to differentiation of HSPC (C). However, the overall number of CD34<sup>+</sup> was highest at Day 7 post-expansion (D). Although the highest frequency of CD34<sup>+</sup>CD90<sup>+</sup>CD45RA<sup>-</sup> cells was observed at Day 5 (E), the highest fold increase in the number of CD34<sup>+</sup>CD90<sup>+</sup>CD45RA<sup>-</sup> cells was observed at Day 7 (F).

Figure 4. Expansion of CD34<sup>+</sup> Umbilical Cord Blood with Our HSPC Expansion Medium at Day 7 post-expansion

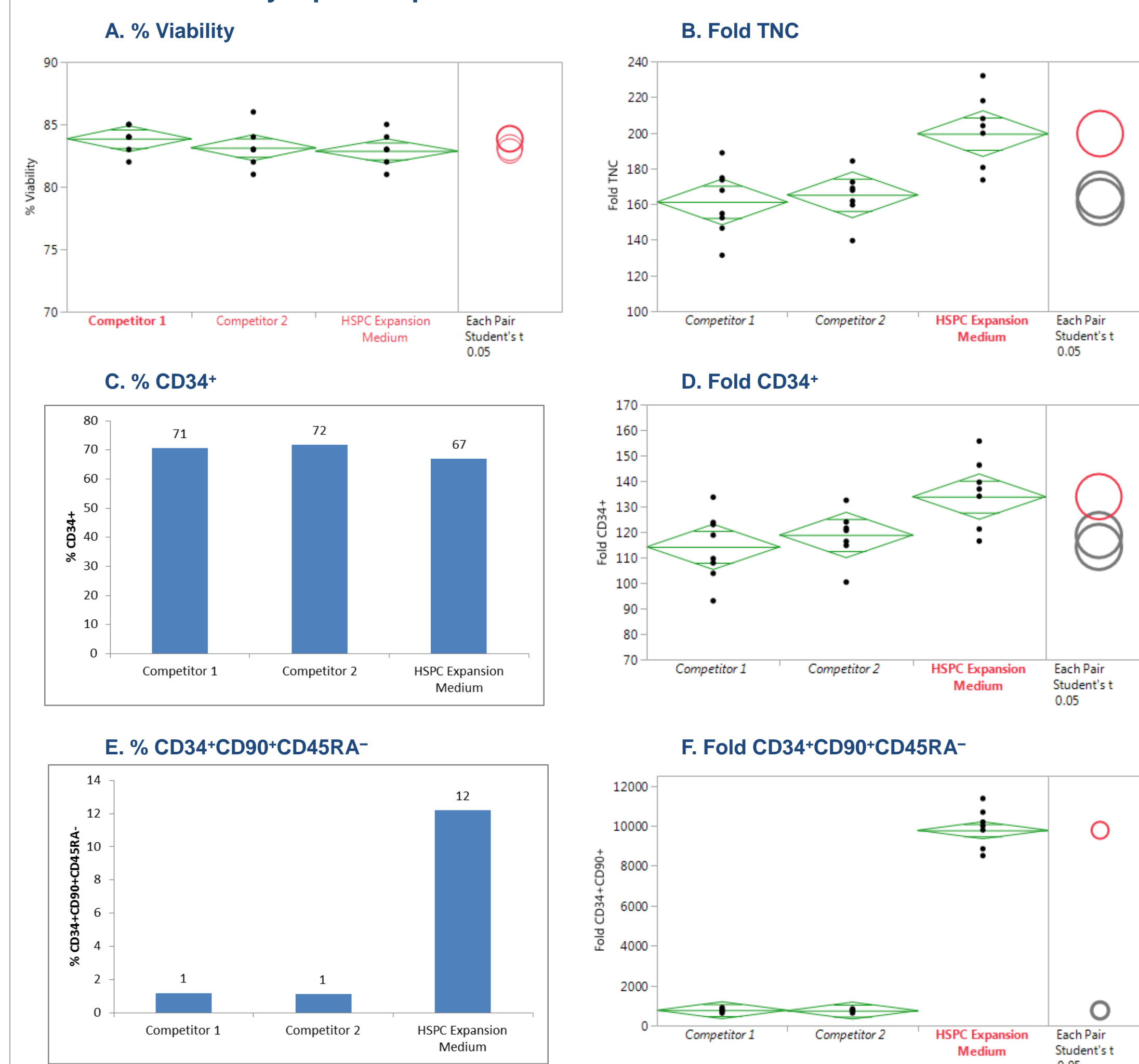


Figure 4: Representative performance of HSPC Expansion Medium when cultured with CD34<sup>+</sup> UCB lot of cells. 7 day expansion with our HSPC Expansion Medium resulted in high % viability, with the highest fold TNC, highest fold CD34<sup>+</sup>, and more importantly significantly high fold expansion of long-term CD34<sup>+</sup>CD90<sup>+</sup>CD45RA<sup>-</sup> compared to available solutions.

Figure 5. Aldehyde dehydrogenase (ALDH) Assay to assess whether expanded cells are bona-fide stem cells

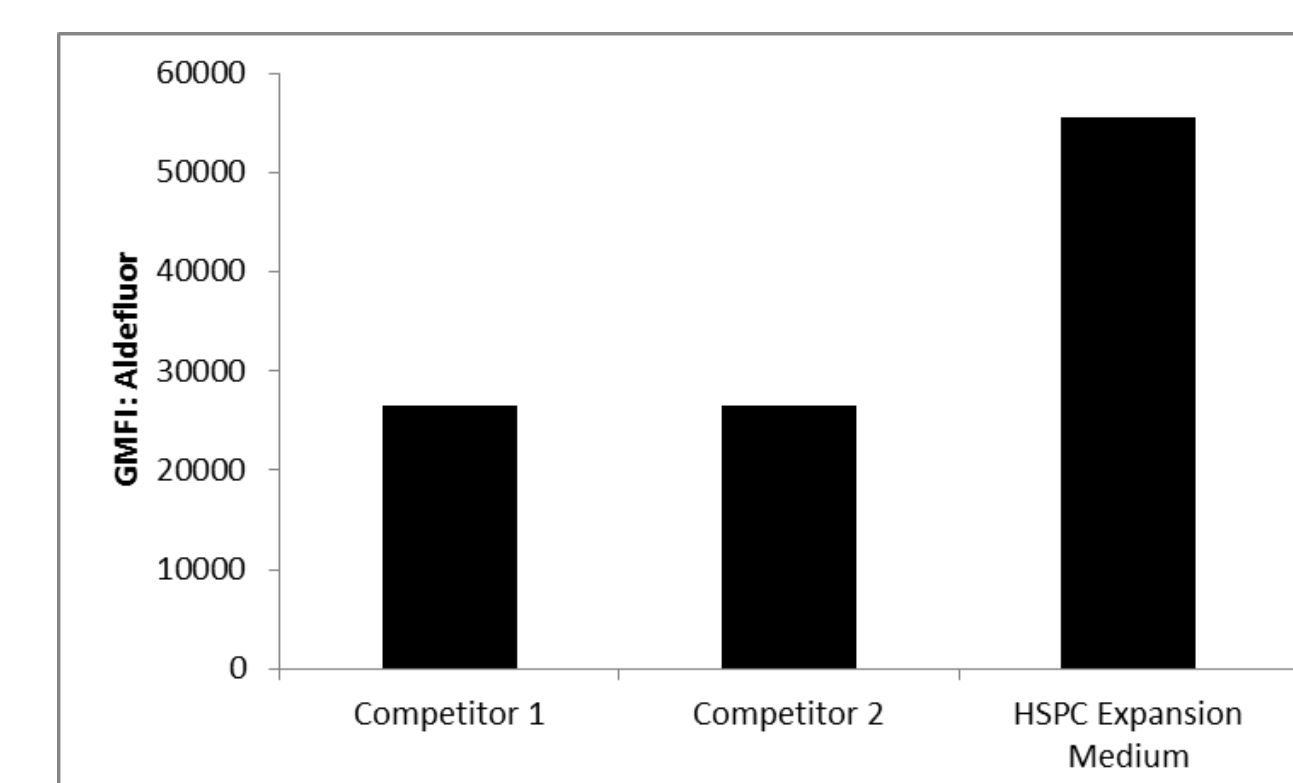


Figure 5: Our HSPC Expansion Medium had the highest ALDH expression as demonstrated by high ALDFLUOR™ staining. This shows that HSPC expanded with our media system contain highest frequency of bona-fide stem cells.

## CONCLUSIONS

- We successfully developed an HSPC Expansion Medium that expands both short term CD34<sup>+</sup> cells, and primitive long-term engraftable hematopoietic stem cells (CD34<sup>+</sup>CD90<sup>+</sup>CD45RA<sup>-</sup>).

## REFERENCES

- Doulatov S, Notta F, Laurenti E, Dick JE. Hematopoiesis: a human perspective. *Cell Stem Cell*, 2012 Feb 3;10(2):120-36.
- Notta F, Doulatov S, Laurenti E, Poeppl A, Jurisica I, Dick JE. Isolation of single human hematopoietic stem cells capable of long-term multilineage engraftment. *Science*, 2011 Jul 8;333(6039):218-21.
- Bhattacharya D, Ehrlich RI, Weissman IL. *European Journal of Immunology*, 2008 Aug; 38(8): 2060–2067.

## ACKNOWLEDGEMENTS

We would like to thank Tracee Crossett, Anne Chen, Melissa Cross, Nicole McCarthy, and Konstantin Slavashevich for their assistance and helpful discussions.

## TRADEMARKS/LICENSING

© 2018 Thermo Fisher Scientific Corporation. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. For Research Use Only. Not for use in diagnostic procedures.