

Stroma-free, serum-free expansion and differentiation of hematopoietic stem and progenitor cells to the T cell lineage

Nooshin Tabatabaei-Zavareh¹, Tim A. Le Fevre¹, Alexander J.Y. Man¹, Evan A. Karas¹, Stephen J. Szilvassy¹, Terry E. Thomas¹, Allen C. Eaves^{1,2}, and Albertus W. Wognum¹

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

Introduction

The use of T cells for cancer immunotherapy and other therapeutic applications relies on the isolation of T cells from peripheral blood and their subsequent activation and expansion in culture. T cells can also be generated from hematopoietic stem and/or progenitor cells (HSPCs) in cord blood (CB) or bone marrow (BM). This approach not only offers a renewable source of T cells, but also provides a model system to study disease mechanisms or validate new drugs that affect T cell development and/or function.

Differentiation of HSPCs to T cells typically requires co-culture with stromal cell lines that have been engineered to express a Notch-ligand. In such cultures, CD34⁺CD38^{-/lo} HSPCs develop into CD7⁺CD5⁺ pro-T cells that further differentiate to T lineage-committed progenitor cells (pre-T cells) characterized by the expression of CD1a. CD7⁺CD1a⁺ pre-T cells can then differentiate to express CD4 and become CD4 immature single-positive (CD4ISP) cells. CD4ISP cells give rise to CD4⁺CD8⁺ double positive (DP) cells. These finally mature into CD4 and CD8 single-positive (SP) CD3⁺TCRαβ⁺ T cells.

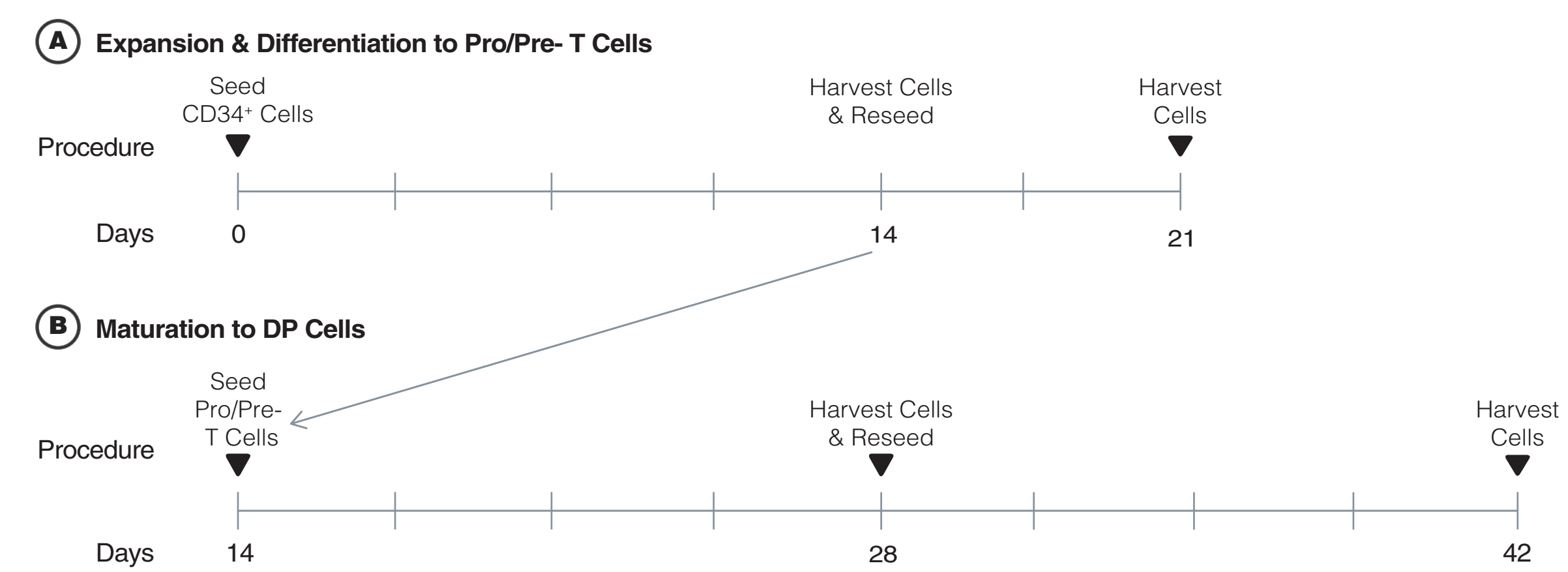
Here we describe a serum-free culture method that recapitulates these differentiation steps in the absence of stromal cells and generates large numbers of functional T lineage cells from a limited number of purified CD34⁺ CB cells. In this system, CD7⁺CD5⁺ pro-T cells were generated with a frequency of 67% (57 - 77%) and a yield of 144 (99 - 188) cells per initial CD34⁺ cell at day 14 (mean & 95% CI; n=25). The frequency and yield of these cells increased to 84% (80 - 88%) and 2130 (1500 - 2761), respectively, when cultured under the same conditions for another week. Pro-T cells were able to differentiate into more mature DP and CD3⁺TCR⁺ T cells when cultured for an additional 4 weeks under different stroma-free culture conditions that support T cell maturation.

Methods

Culture Protocol

CD34⁺ cells were enriched from human CB samples by depleting mature cells using RosetteSep™ followed by EasySep™ CD34⁺ positive selection. The isolated CD34⁺ cells were plated at 1x10⁴ cells/mL in StemSpan™ SFEM II medium supplemented with T Cell Progenitor Expansion Supplement (containing SCF, TPO, Flt3L, and IL-7) onto plates pre-coated with StemSpan™ T Cell Progenitor Differentiation Coating Material (StemSpan™ T Cell Progenitor Differentiation Kit). Every 3 - 4 days a half medium exchange was performed. The cells were cultured for 14 days, after which they were harvested, counted and re-plated at 1x10⁵ cells/mL onto freshly coated plates for an additional 7 days of culture. T lineage cells were harvested on day 14 or 21 for analysis and/or further maturation. This protocol is detailed in Figure 1A. For further maturation the pro/pre-T cells harvested on day 14 were seeded at 1x10⁵ cells/mL on freshly coated plates with the same attachment substrate in a modified T Cell Progenitor Maturation medium containing IL-7 and Flt3L. After 14 days, cells were harvested, counted and re-plated at 2 x 10⁵ cells/mL onto freshly coated plates for an additional 14 days of culture. Matured DP T cells were harvested after a total of 6 weeks of culture (Figure 1B).

FIGURE 1: Workflow and Culture Protocol

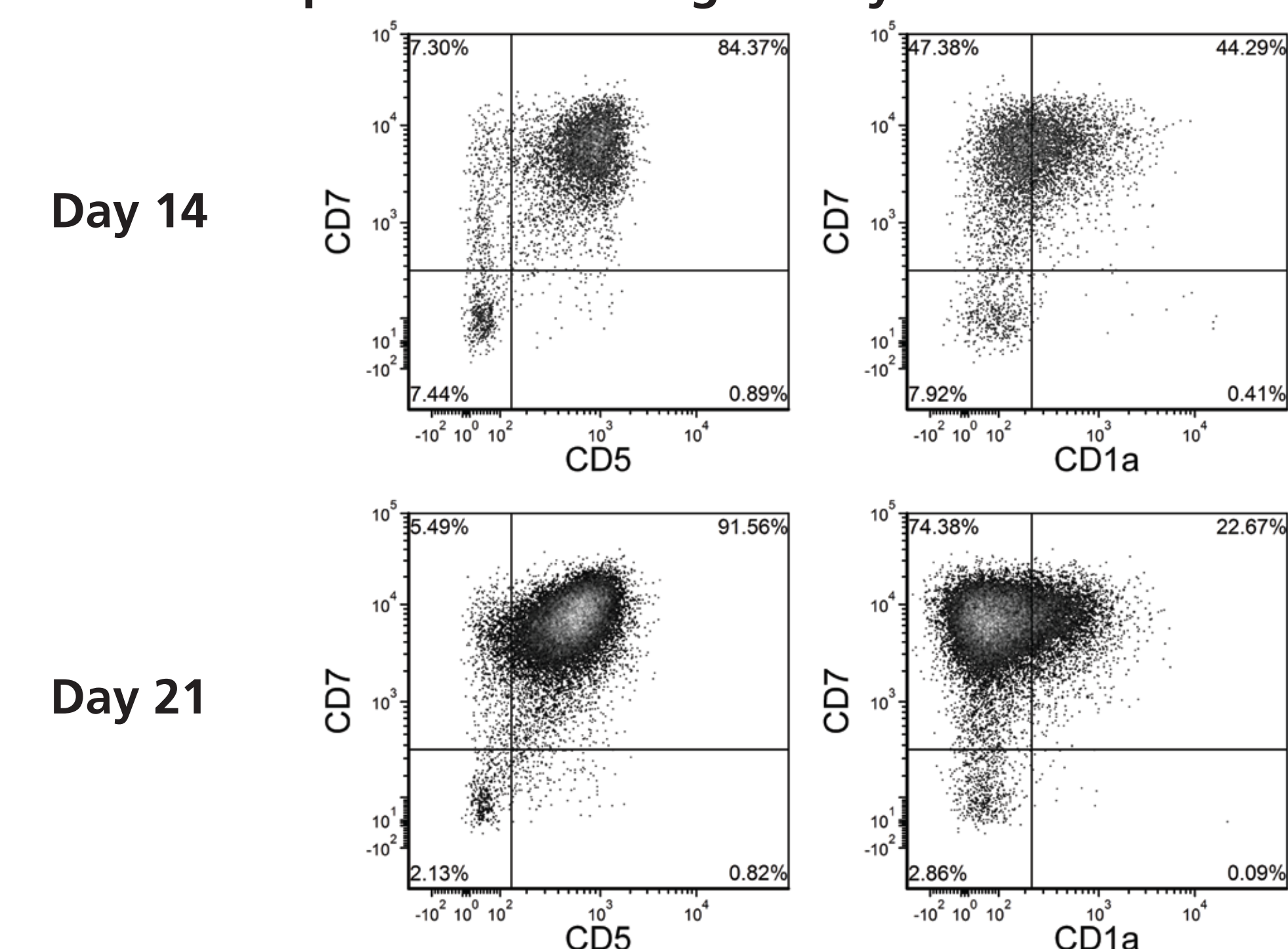


Assessment of T Lineage Cells

Harvested cells were counted and analyzed by flow cytometry for the expression of T lineage markers including CD7, CD5, CD1a, CD4, CD8, CD3, TCRαβ and TCRγδ. Dead cells were excluded by light scatter profile and 7-AAD staining. The number of pro-T (CD5⁺CD7⁺), pre-T (CD7⁺CD1a⁺), CD4ISP (CD4⁺CD3⁻TCR⁻), DP (CD4⁺CD8⁺) and CD3⁺TCRαβ⁺ T cells (Figure 3 and 4) was calculated based on the fraction of cells counted expressing the specified markers.

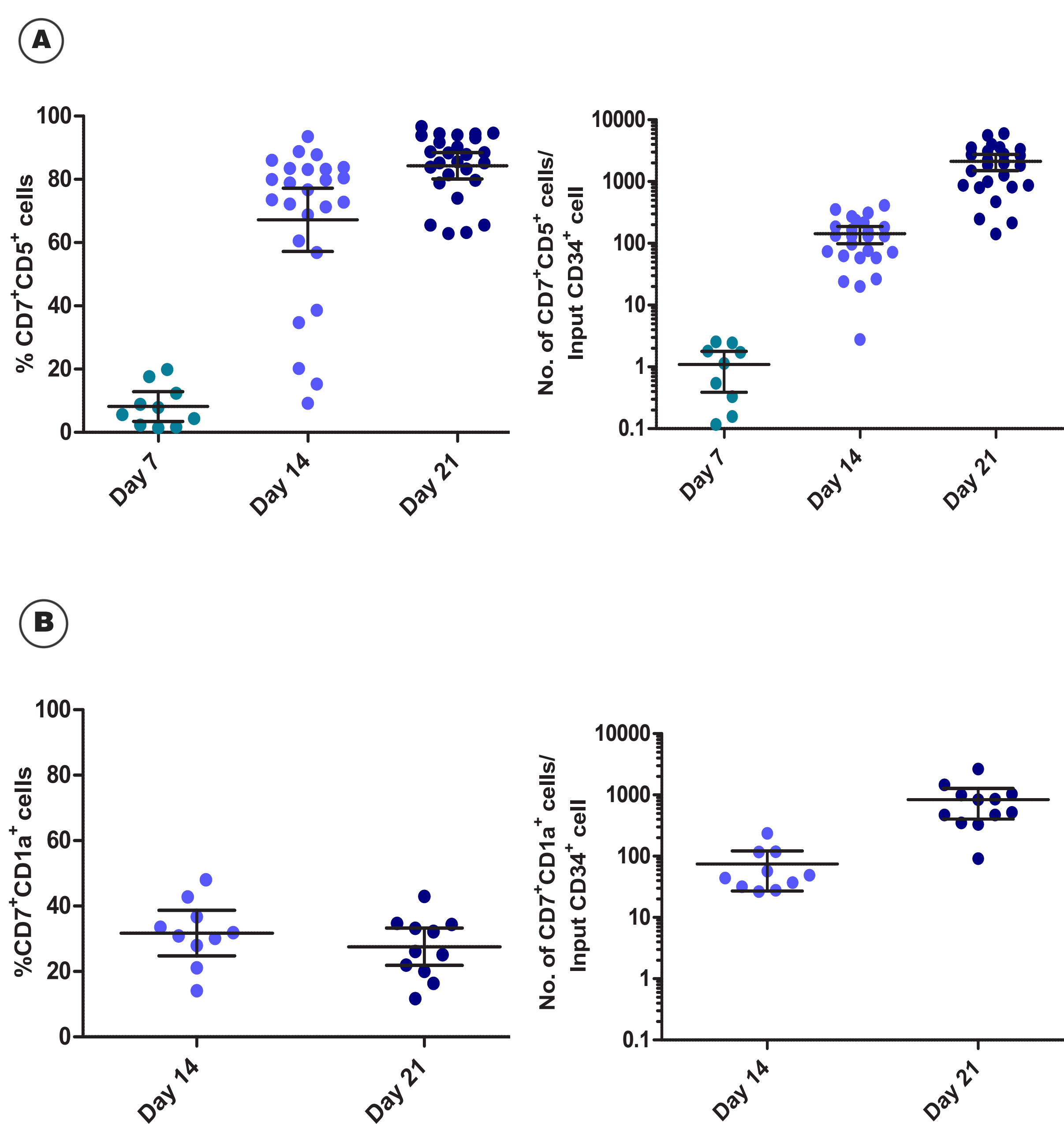
Results

FIGURE 2: CD34⁺ CB cells differentiate into CD7⁺CD5⁺ pro-T and CD7⁺CD1a⁺ pre-T cells during 21 days of culture



CD34⁺ CB cells (1,000 cells/well) were cultured for 3 weeks. Cells were analyzed by flow cytometry for the expression of CD7, CD5 and CD1a on days 14 and 21.

FIGURE 3: StemSpan™ T Cell Progenitor Differentiation Kit promotes expansion of CD34⁺ CB cells and their differentiation into CD7⁺CD5⁺ pro-T and CD7⁺CD1a⁺ pre-T cells

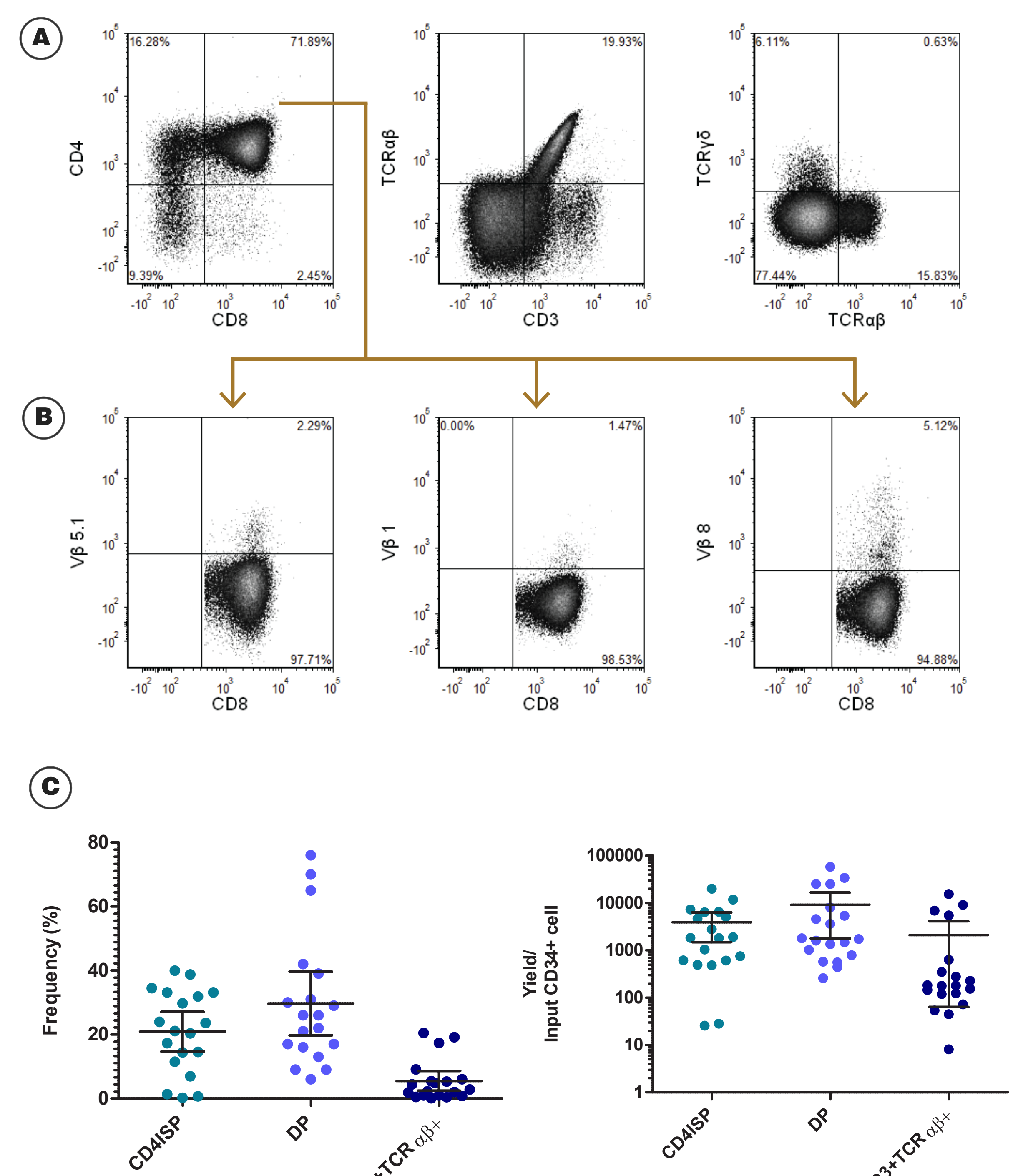


CD34⁺ CB cells were cultured for the indicated times as shown in Figure 1A, and then counted and analysed by flow cytometry as described in Methods. The percentage and number of (A) CD7⁺CD5⁺ cells and (B) CD7⁺CD1a⁺ cells generated are shown for 10 - 26 independent experiments. Horizontal lines indicate the mean. Vertical lines indicate 95% confidence interval. (A) The frequency of CD7⁺CD5⁺ pro-T cells increased each week up to, on average, 84% on day 21. The number of pro-T cells also increased ~10 - 100 fold every week, resulting in an average of ~2000 pro-T cells per initial CD34⁺ cell on day 21. (B) On day 21 on average 28% of the cells express CD7 and CD1a, indicating further differentiation into pre-T cells. The yield of CD7⁺CD1a⁺ cells on day 21 was ~800 per initial CD34⁺ cell.

Conclusions

- CD34⁺ HSPCs from CB proliferate and differentiate efficiently into T cell progenitors in the absence of serum and stromal cells when cultured for 2 - 3 weeks using serum-free StemSpan medium and T Cell Progenitor Differentiation Supplements.
- This culture system can generate >2000 CD7⁺CD5⁺ pro-T cells with ~80% frequency (and ~800 pre-T cells) per original CD34⁺ CB cell during 3 weeks of culture.
- Pro- and pre-T cells can differentiate further into DP and CD3⁺TCR⁺ T cells during 4 weeks of continued stroma-free culture using a T Cell Progenitor Maturation Supplement.
- The overall yield of CD3⁺TCRαβ⁺ T cells is ~2000 per initial CD34⁺ CB cell (with average frequency of 6%) after a total of 42 days of culture.

FIGURE 4: Pro-T cells differentiate into CD4ISP, DP and CD3⁺TCR⁺ T cells during continued stroma-free culture



50,000 pro/pre-T cells generated after 14 days of culture as described in Figure 1A were replated in StemSpan™ SFEM II medium supplemented with a T Cell Progenitor Maturation Supplement (containing Flt3L and IL-7) in 24-well plates coated with StemSpan™ T Cell Differentiation Coating Material and cultured for another 4 weeks (Figure 1B). (A & B) Cells were then analyzed by flow cytometry for the expression of CD3, CD4, CD8, TCRαβ and TCRγδ. (B) Cells were also stained with antibodies that recognize variants of TCR Vβ chain to assess polyclonality of TCR. These cells were first gated on DP cells. (C) Cell counts were also obtained. Data shows the mean with 95% CI of 19 experiments. On average 21% CD4⁺CD8⁻ CD4ISP, 30% CD4⁺CD8⁺ DP and 6% CD3⁺TCRαβ⁺ cells arise in these stroma-free maturation cultures with a yield of ~4000 CD4ISP, 9000 DP and 2000 CD3⁺TCRαβ⁺ cells, respectively, per original CD34⁺ cell. The majority of CD3⁺TCRαβ⁺ cells are DP.