

# Improved *Ex Vivo* Expansion of Hematopoietic Progenitor Cells and their Differentiation into Megakaryocytes and Erythroid Cells in a Novel Serum-Free Medium, StemSpan SFEM II

Bert Wognum<sup>1</sup>, Donna DeGeer<sup>1</sup>, Ning Yuan<sup>1</sup>, Terry Thomas<sup>1</sup> & Allen Eaves<sup>1,2</sup>

<sup>1</sup>STEMCELL Technologies Inc, Vancouver, BC, Canada

<sup>2</sup>Terry Fox Laboratory, Vancouver, BC, Canada

## Abstract

We developed a new serum-free medium, StemSpan SFEM II, for *ex vivo* expansion and lineage-specific differentiation of human hematopoietic progenitor cells (HPCs). SFEM II is an enhanced version of the original SFEM medium, which has been supplemented with a chemically-defined lipid formulation. To compare the performance of SFEM II to other media formulated for HPC expansion, cord blood (CB) CD34<sup>+</sup> cells (10<sup>4</sup>/mL) were cultured in SFEM II and, as controls, in SFEM I, H3000, and animal component-free (ACF) medium. The cultures were supplemented with Flt-3 Ligand (FL), stem cell factor (SCF), interleukin (IL)-3, and IL-6. After 7 days, cells were counted, analyzed for CD34 expression and in some experiments assayed to measure colony-forming units (CFUs). Cell viability was high in all cultures (range 93 to 99%; n=10). Total nucleated cell (TNC) expansion in H3000, SFEM I, and ACF media averaged 40-fold (range 14 to 76-fold) and was not significantly different between the three media (P>0.1; paired t-test). TNC expansion in SFEM II medium was significantly higher (on average 60-fold expansion; range 23 to 101-fold; p<0.01; n=10). CD34<sup>+</sup> cell output was also highest in SFEM II, but only the differences with H3000 and SFEM I were significant (p<0.01). CFC expansion showed similar differences, with highest expansion of BFU-E, CFU-GM, and CFU-GEMM numbers in SFEM II, and lowest in H3000. The capacity of SFEM II and SFEM I to support megakaryocytic (Mk) and erythroid progenitor expansion and differentiation of CD34<sup>+</sup> cells was tested in cultures that were refed on days 4, 7, and/or 10 to maximize expansion and prevent overgrowth, and analysed on day 14. Using a Mk-selective cytokine cocktail [thrombopoietin (TPO), SCF, IL-6, and IL-9], expansion of CD45<sup>+</sup>CD41<sup>+</sup> cells ranged from 13 to 124-fold in eight different CB samples cultured in SFEM I and from 18 to 326-fold in SFEM II, with on average 2-fold higher expansion in SFEM II (p<0.01; n=8). In erythroid-selective cytokines (SCF, IL-3, erythropoietin) and with a glucocorticoid receptor agonist (dexamethasone or hydrocortisone), both media strongly supported production of erythroid cells with 200 to 6,500-fold and 900 to 7,000-fold expansion of cells expressing glycophorin-A (GpA) and/or CD71 in SFEM I and SFEM II, respectively (p<0.05; n=7). These data indicate that StemSpan SFEM II is a superior medium to support the expansion of CB CD34<sup>+</sup> progenitors and to promote their differentiation into the Mk and erythroid lineages. SFEM II should prove useful for the continued development of new cellular therapy applications of culture-expanded HSPCs and mature blood cells.

## Materials and Methods

- Cord blood CD34<sup>+</sup> cells isolated by EasySep HPC Enrichment Kit (STEMCELL, Cat. # 19356) or EasySep Human Cord Blood CD34<sup>+</sup> Selection Kit (STEMCELL, Cat. # 18096) were plated at 10,000 CD34<sup>+</sup> cells per mL.
- Culture conditions for TNC and CD34<sup>+</sup> cell expansion: FL, SCF, IL-3, and IL-6 (CC100 cytokine cocktail); 7 days of culture.
- Culture conditions for Mk expansion and differentiation: TPO, SCF, IL-6, and IL-9 (CC220 cytokine cocktail); 14 days of culture with addition of fresh medium on day 7.
- Culture conditions for erythroid expansion and differentiation: SCF, IL-3, and EPO and either 1 μM dexamethasone or hydrocortisone; 14 days of culture with medium changes or replating on days 4, 7, and 10.
- Analysis after culture: Total and viable cell counting and immunophenotyping by flow cytometry; CFU assay by replating in semisolid medium (MethoCult H4034).

## Results

*Total and CD34<sup>+</sup> cell expansion in 14-day cultures in four StemSpan media supplemented with FL, SCF, IL-3, and IL-6 (CC100 cytokine cocktail).*

Approximately 10-11% of cells in SFEM I, SFEM II, and ACF media retained CD34<sup>+</sup> expression, while the %CD34<sup>+</sup> cells in StemSpan H3000 was 2-fold lower (5.6%; p<0.01, n=10; **Figure 1**). TNC expansion in H3000, SFEM I, and ACF averaged 40-fold (range 14 to 76-fold; p>0.1, paired t-test) and was significantly higher in SFEM II (average 60-fold; range 23 to 101-fold; p<0.01; n=10) (**Figure 2A**). CD34<sup>+</sup> cell expansions in SFEM II and ACF were similar but higher than in SFEM I and H3000 (p<0.01; **Figure 2B**). Expansion of CFUs showed similar differences and similar results were obtained in cultures of bone marrow CD34<sup>+</sup> cells and after stimulation with SCF, FLT3-L, and TPO (CC110 cytokine cocktail; data not shown).

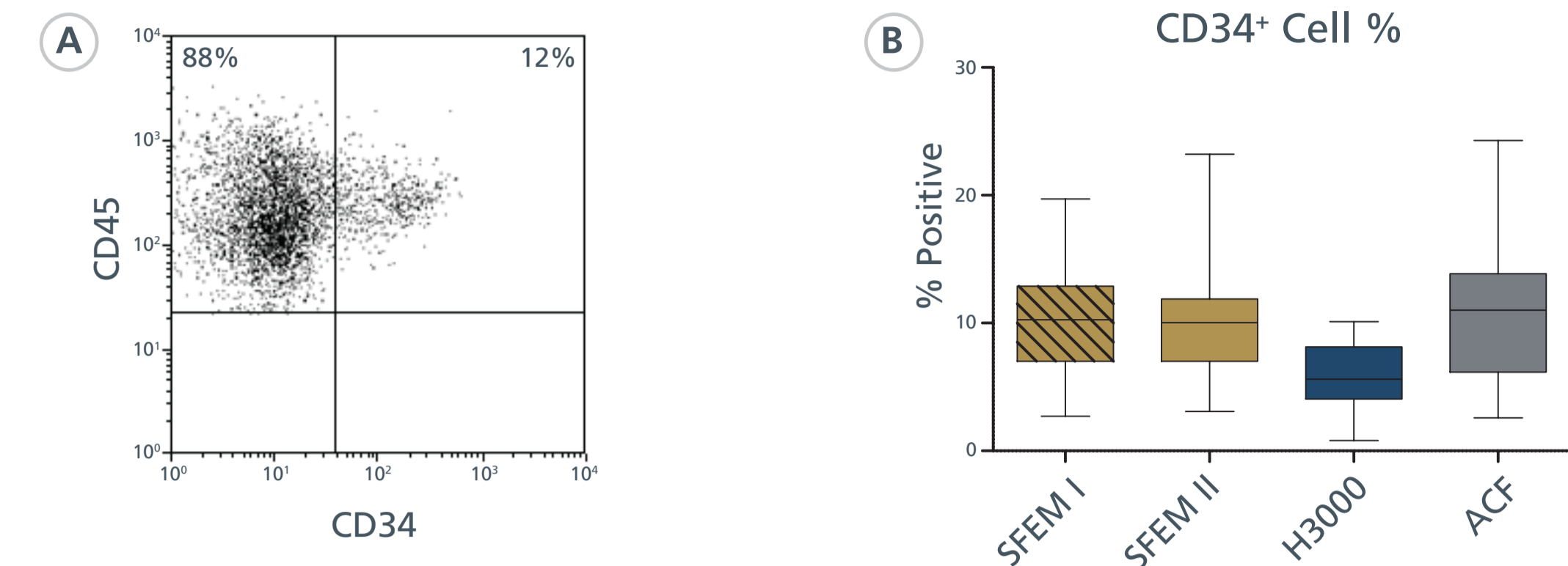
*Megakaryocyte cultures in SFEM I and SFEM II supplemented with TPO, SCF, IL-6, and IL-9 (CC220 cytokine cocktail).*

Both media yielded Mk lineage cells with similarly high purity, i.e., %CD45<sup>+</sup>CD41a<sup>+</sup> cells = 90 ± 6 and 88 ± 6 in SFEM I and SFEM II, respectively (p=0.2; n=8; **Figure 3A,B**). Average Mk cell output in SFEM II (98 CD45<sup>+</sup>CD41a<sup>+</sup> cells per input CD34<sup>+</sup> cell) was two-fold higher than in SFEM I (48 Mk per input CD34<sup>+</sup> cell; p<0.01; n=8; **Figure 3C**). Mk expansion in SFEM I was promoted by supplementation with low-density lipoproteins (LDL, 20 μg/mL), but Mk expansion in SFEM II was 1.6 ± 0.2 fold higher than in SFEM I + LDL (range 1.3 to 1.8-fold; p<0.05; n=4; **Table 1**). Approximately 90% of the CD41a<sup>+</sup> cells expressed the late Mk marker CD42. CD41a<sup>+</sup>CD42<sup>+</sup> platelet-like particles with similar light scatter properties as fresh blood platelets were also detectable, indicating that the media supported terminal Mk differentiation (data not shown).

*Erythroid cultures in SFEM I and SFEM II supplemented with SCF, IL-3, EPO, and glucocorticoid receptor agonist (dexamethasone or hydrocortisone at 1 μM).*

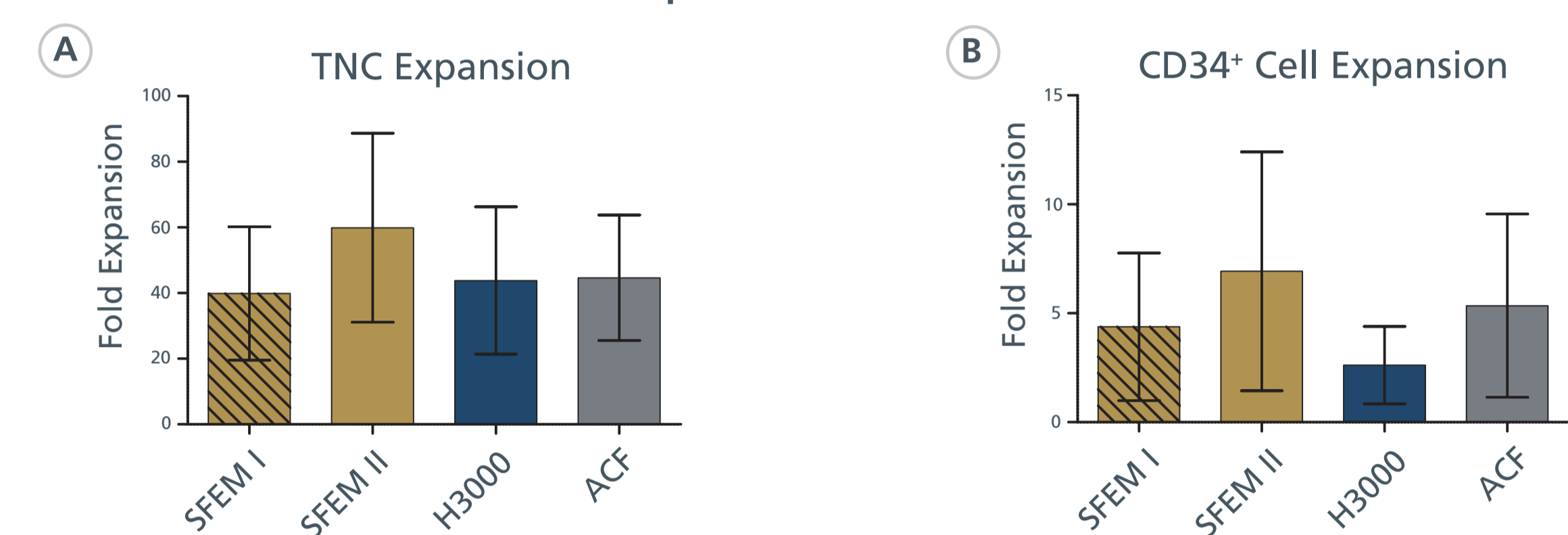
Both media yielded similar erythroid cell frequencies: average ±SD = 76 ± 13% and 81 ± 6% in SFEM I and II, respectively (p=0.3; n=7; **Figure 4A,B**). Approximately 40-60% of cells were CD71<sup>+</sup>GpA<sup>+</sup> erythroblasts, while CD71<sup>+</sup>GpA<sup>-</sup> pro-erythroblasts and CD71<sup>-</sup>GpA<sup>+</sup> normoblasts were much less frequent (**Figure 4B**). Cell output in SFEM II (average: 2,800 erythroblasts per input CD34<sup>+</sup> cell; range 900 to 7,000) was ~20% higher than in SFEM I (average: 2,300; range 200 to 6,500; p<0.05; n=7; **Figure 4C**). Cumulative cell output in three cultures that were maintained for 26 days ranged between 10<sup>5</sup> and 10<sup>6</sup> erythroblasts per input CD34<sup>+</sup> cell (data not shown).

**FIGURE 1: CD34 expression after 7 days of culture of purified CB CD34<sup>+</sup> cells in four different StemSpan media**



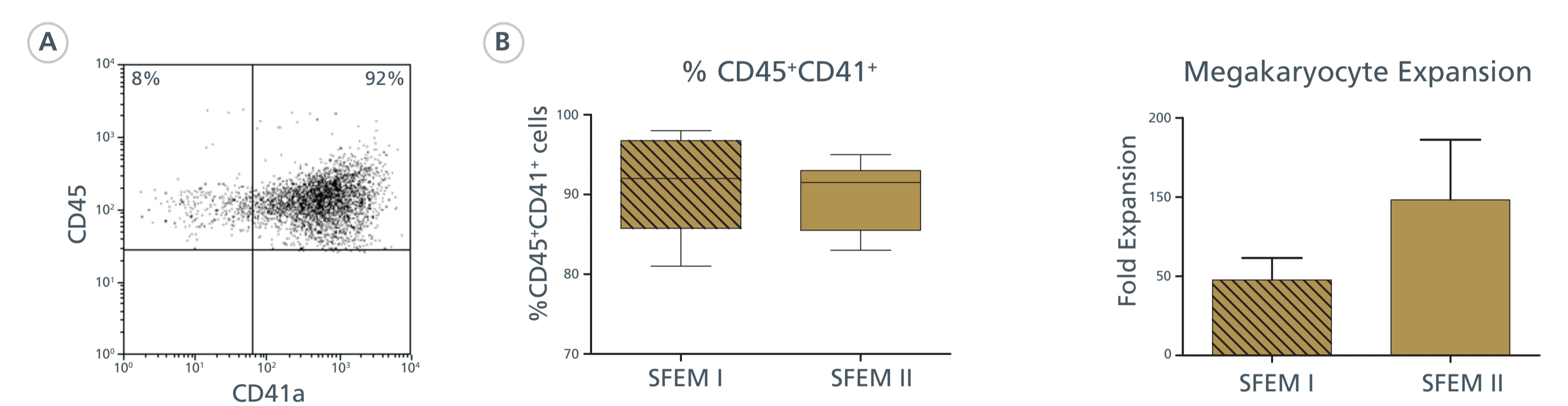
**A**) Flow cytometry dot plot showing CD45 and CD34 expression after 7 days of culture of CB CD34<sup>+</sup> cells in medium supplemented with FL, SCF, IL-3, and IL-6 (CC100 cytokine cocktail). **B**) Percentage of CD34<sup>+</sup> cells as measured by flow cytometry as depicted in the example plot shown in A. Culture of CB CD34<sup>+</sup> cells in SFEM I, SFEM II, and ACF media resulted in similar frequencies of CD34<sup>+</sup> cells (p>0.5; paired t-test; n=10), which were ~2-fold higher than in H3000 medium (p<0.01). Box and whisker plot show median, 25-75 percentile and full data range.

**FIGURE 2: Total and CD34<sup>+</sup> cell expansion after 7 days of culture of purified CB CD34<sup>+</sup> cells in four different StemSpan media**



Average fold increase of **A**) Total nucleated cells (TNC) and **B**) CD34<sup>+</sup> cells after 7 days of culture. Bars represent standard deviation (n=10). SFEM II supported, on average, ~50% higher TNC expansion compared to the other three media (p<0.01). SFEM II and ACF media supported similar ~8-fold CD34<sup>+</sup> cell expansion (p>0.5; paired t-test; n=10), which was ~2-fold higher than in H3000 medium (p<0.01).

**FIGURE 3: Megakaryocyte outgrowth of CB CD34<sup>+</sup> cells in two StemSpan SFEM media**

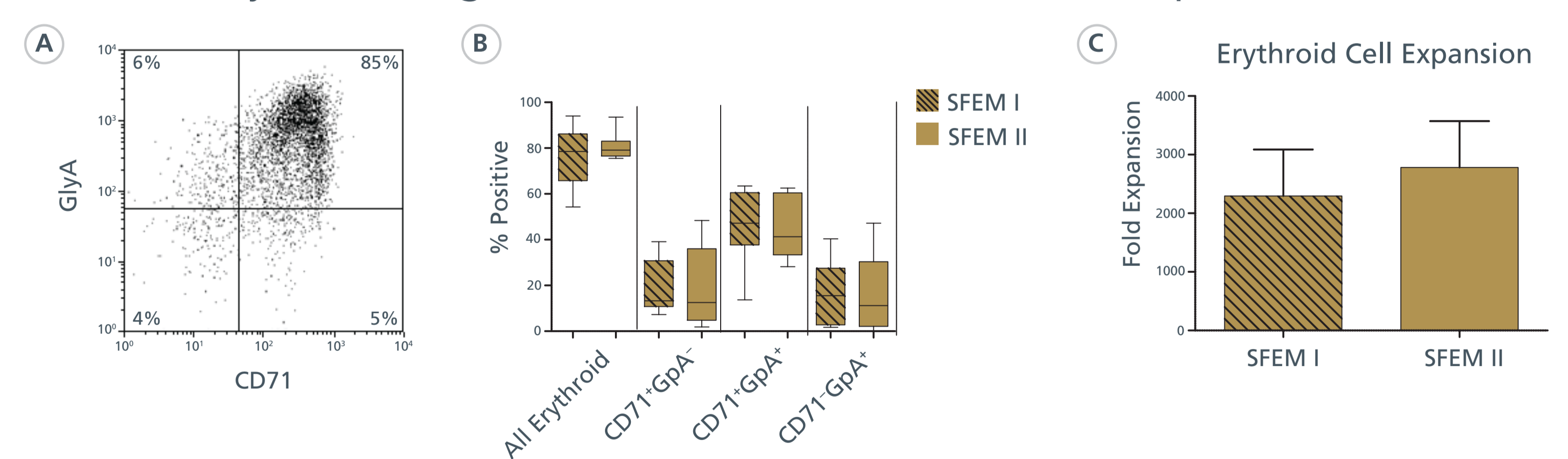


**A**) Example of CD45 and CD41 expression after 14 days of culture of CB CD34<sup>+</sup> cells under Mk-specific conditions (TPO, SCF, IL-6, and IL-9, CC220 cytokine cocktail). **B**) Percentage of CD45<sup>+</sup>CD41<sup>+</sup> Mk was measured by flow cytometry as depicted in the sample plot shown in A. **C**) Average fold Mk expansion, expressed as number of CD45<sup>+</sup>CD41<sup>+</sup> cells generated per input CD34<sup>+</sup> cell. Bars represent standard deviation (n=8).

**TABLE 1: SFEM II supports higher Mk output in cultures of CB CD34<sup>+</sup> cells compared to SFEM I supplemented with low-density lipoproteins**

Exp	Fold expansion of CD45 <sup>+</sup> CD41 <sup>+</sup> cells		
	SFEM I	SFEM I + LDL	SFEM II
1	16.0	19.7	35.7
2	20.7	25.8	44.3
3	47.0	53.2	67.6
4	124.1	213.6	326.4
Average	52.0	78.1	118.5
SD	50.0	91.5	139.2

**FIGURE 4: Erythroid outgrowth of CB CD34<sup>+</sup> cells in two StemSpan SFEM media**



**A**) Example of CD71 and GpA expression after 14 days of culture of CB CD34<sup>+</sup> cells under erythroid-specific culture conditions (SCF, IL-3, EPO, dexamethasone). **B**) Percentage of all erythroid cells and of erythroid subsets identified on the basis of CD71 and GpA expression as shown in A. **C**) Average fold expansion of all erythroid cells (CD71<sup>+</sup> and/or GpA<sup>+</sup>) expressed as number of cells generated per input CD34<sup>+</sup> cell. Bars represent standard deviation (n=7).

## Conclusions

- **StemSpan SFEM II supports stronger *ex vivo* expansion of human CB and BM hematopoietic cells, with higher numbers of nucleated cells and CD34<sup>+</sup> cells in short-term (7 days) cytokine supplemented cultures than original StemSpan SFEM (SFEM I), a xeno-free medium (StemSpan H3000) and an animal and human serum protein-free medium (StemSpan-ACF).**
- **StemSpan SFEM II strongly supports megakaryocyte (Mk) expansion and differentiation of CB CD34<sup>+</sup> cells under Mk-specific culture conditions, resulting in ~100-fold cell expansion and generation of ~90% pure Mk lineage cells after 14 days of culture.**
- **StemSpan SFEM II also strongly supports erythroid expansion and differentiation of CB CD34<sup>+</sup> cells under erythroid-specific culture conditions, resulting in large cell expansion (>10<sup>3</sup> and >10<sup>5</sup> fold after 14 and 26 days, respectively) and generation of highly pure populations of CD71<sup>+</sup> and/or Glycophorin-A<sup>+</sup> erythroid lineage cells.**