

# Rapid and Efficient Positive Selection of Hematopoietic Stem and Progenitor Cells from Adult Mouse Bone Marrow

M Fairhurst<sup>1</sup>, SM Woodside<sup>1</sup>, AW Wognum<sup>1</sup>, CL Miller<sup>1</sup>, AC Eaves<sup>2</sup> and TE Thomas<sup>1</sup>

<sup>1</sup>StemCell Technologies Inc, Vancouver, Canada, and <sup>2</sup>Terry Fox Laboratory, BC Cancer Agency, Vancouver, Canada

## Introduction

No antigenic marker has yet been shown to be uniquely expressed by murine hematopoietic stem cells (HSCs) or their immediate progeny. However, most of these cells express Sca-1 (Ly6A/E) and c-kit (CD117), neither of which is expressed on the majority of cells in murine hematopoietic tissues, including adult bone marrow (BM). Sca-1 and c-kit are therefore useful as single markers for obtaining enriched populations of primitive hematopoietic cells from this source. Here we describe a preparative immunomagnetic enrichment procedure that allows the rapid and direct selection of c-kit<sup>+</sup> or Sca-1<sup>+</sup> cells from adult mouse BM using FACS-compatible superparamagnetic nanoparticles. These results demonstrate the utility of this approach for the rapid enrichment of primitive murine hematopoietic cells and show that the functional integrity of repopulating cells *in vivo* is not compromised by the labeling and isolation procedure.

## Methods

**Bone marrow cells.** 10<sup>8</sup> BM cells were used for each protocol. EasySep™ magnetic positive selection was used to select c-kit<sup>+</sup> and Sca-1<sup>+</sup> cells. Lineage-positive cells were removed (lineage-depleted) using a cocktail of antibodies (CD5, CD11b, CD45R, Gr-1, 7-4, Ter119) and SpinSep™ dense particle sedimentation (StemCell Technologies Inc).

**Analysis of enriched cells.** The positively selected cells are labeled with the PE-conjugated antibody during the EasySep™ selection protocol and purity can be assessed by flow cytometry directly or after co-staining with either anti-c-kitAPC or anti-Sca-1Alexa647 and a cocktail of lineage markers [(CD3, CD11b, CD45R, Gr-1, Ter119)FITC].

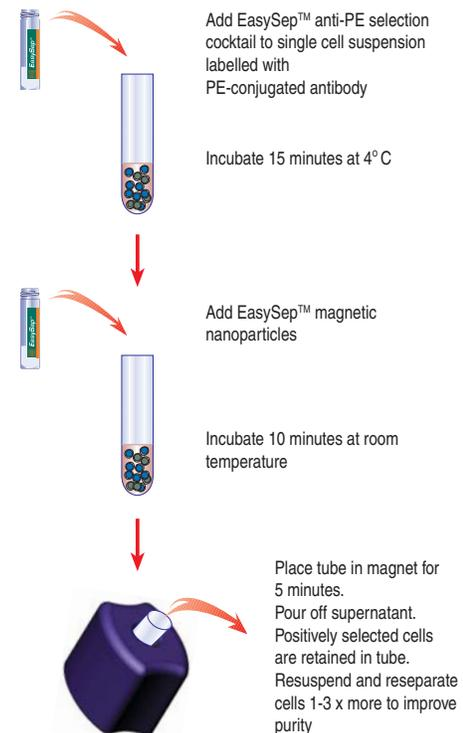
**Colony forming cell (CFC) assay.** Pre- and post-enriched BM cells were plated in MethoCult™ GF 3434. Unseparated cells were plated at 10<sup>4</sup>/dish. Enriched cells were plated at 1000/dish. Colonies were counted on day 10-12. Total CFCs were defined as BFU-E + CFU-GEMM + CFU-GM.

**Competitive repopulating unit (CRU) assay.** C57Bl/6 (Ly5.2<sup>-</sup>) donor cells were injected into groups of irradiated (9 Gy) C57Bl/6/Pep3B (Ly5.1<sup>+</sup>) recipients together with 10<sup>5</sup> C57Bl/6/Pep3B (Ly5.1<sup>+</sup>) BM cells. The following cell doses were used per recipient:

- A: unseparated (15000, 45000, 90000, 135000);
- B: SpinSep™ lineage-depleted (4000, 12000, 24000, 36000);
- C: c-kit<sup>+</sup> 4x separated (1500, 4500, 9000, 13500);
- D: c-kit<sup>+</sup> 2x separated (3000, 9000, 18000, 27000);
- E: Sca-1<sup>+</sup> (2500, 7500, 15000, 22500).

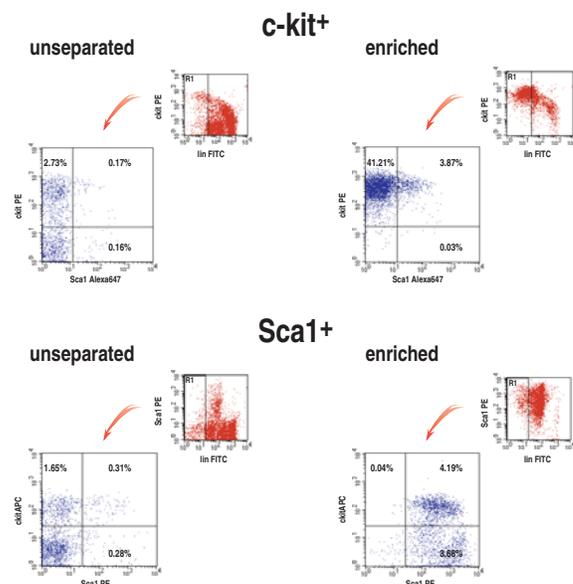
The assay was terminated at 4 months and the presence of donor-derived (Ly5.2<sup>-</sup>) myeloid (Gr1<sup>+</sup>) and lymphoid (CD3<sup>+</sup>/CD19<sup>+</sup>) WBCs was assessed (reference 1).

Figure 1. EasySep™ procedure for cell selection



## Results

Figure 2. Typical FACS plots of unseparated and c-kit<sup>+</sup> or Sca-1<sup>+</sup> mouse BM cells after EasySep™ immunomagnetic positive selection



Plots are gated on PI<sup>-</sup>, lineage [(CD3, CD11b, CD45R, Gr-1, Ter119) FITC] negative cells = R1



StemCell Technologies  
Tel: 604.877.0713 • Fax: 604.877.0704  
Toll Free Tel: 1.800.667.0322 • Toll Free Fax: 1.800.567.2899  
E-mail: info@stemcell.com  
www.stemcell.com

**Table 1. Purity and recovery of c-kit<sup>+</sup> or Sca-1<sup>+</sup> mouse BM cells after EasySep™ immunomagnetic positive selection**

target cell	phenotype	n	% purity start	% purity enriched	% recovery	fold enrichment	n	% purity enriched	% recovery	fold enrichment
				2x separated*				4x separated*		
c-kit <sup>+</sup>	c-kit <sup>+</sup>	6	20±1	68±4	34±7	3.7±0.3	12	92±1	12±2	4.9±0.4
	lin-c-kit <sup>+</sup> Sca-1 <sup>+</sup>	3	0.17±0.03	1.5±0.2	58±7	9.2±0.9	3	3.9±0.6	43±5	24±5
Sca-1 <sup>+</sup>	Sca-1 <sup>+</sup>	7	10.8±0.7	70±2	58±4	6.7±0.4	13	94±1	30±2	9±1
	lin-c-kit <sup>+</sup> Sca-1 <sup>+</sup>	6	0.43±0.04	3.6±0.4	69±5	7.6±0.8	9	4.3±0.5	37±5	9.9±0.5

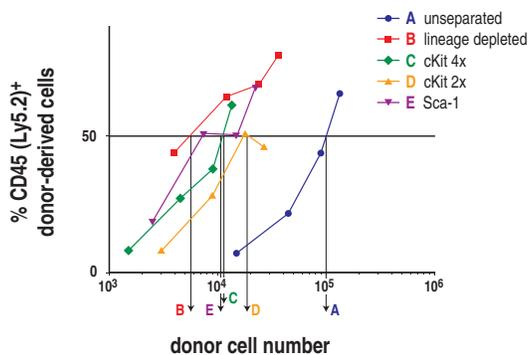
values are means ± 1 sem; n = number of experiments  
\*2 or 4 rounds of separation

**Table 2. Recovery and fold enrichment of CFCs from mouse BM enriched for c-kit<sup>+</sup> or Sca-1<sup>+</sup> cells using EasySep™ immunomagnetic positive selection**

target cell	n	CFC per 1000 cells	total CFC % recovery	fold enrichment	n	CFC per 1000 cells	total CFC % recovery	fold enrichment	
unseparated	12	3.4±0.3			12	3.4±0.3			
			2x separated*				4x separated*		
c-kit <sup>+</sup>	6	32±6	77±6	9±2	11	70±5	57±5	26±4	
Sca-1 <sup>+</sup>	6	23±3	54±7	6±1	12	27±3	35±4	10±2	

Values are means ± 1 sem; n = number of experiments  
\*2 or 4 rounds of separation

**Figure 3. Percent repopulation in irradiated recipients given unseparated or enriched cell fractions**



Engrafted mice showed lymphoid and myeloid repopulation. Data points represent the average % donor-derived repopulation for 3 animals. Number of donor cells required for 50% repopulation (A, B, C, D, E).

**Table 3. Frequency, fold enrichment and recovery of CRU from unseparated, lineage depleted, c-kit<sup>+</sup> or Sca-1<sup>+</sup> selected mouse BM cells**

cells	number of donor cells at 50% engraftment level*	CRU frequency**	fold enrichment***	% recovery CRU****
A: unseparated	100,000	1/16,000	-	-
B: lineage depleted	5,400	1/864	19	13
D: c-kit <sup>+</sup> 2x separated	18,000	1/2,880	6	74
C: c-kit <sup>+</sup> 4x separated	11,000	1/1,760	9	29
E: Sca-1 <sup>+</sup>	10,500	1/1,680	10	59

Results of one experiment with 3 mice per group. c-kit<sup>+</sup> and Sca-1<sup>+</sup> cells were selected using EasySep™ immunomagnetic positive selection. Lineage-depleted cells obtained using SpinSep™ dense particle sedimentation.  
\* from Fig. 3: A, B, C, D and E  
\*\*CRU frequency = 1/(number donor-derived cells at 50% repopulation level per 6.25 CRU)  
(assuming that 8% repopulation = 1 CRU; therefore, 50% repopulation = 6.25 CRU)  
\*\*\* (CRU frequency in enriched fraction)/(CRU frequency in unseparated fraction)  
\*\*\*\* [(total number of cells x CRU frequency)<sub>post separation</sub> / (total number of cells x CRU frequency)<sub>unseparated</sub>] x 100

## Conclusions

- Hematopoietic stem cells and progenitor cells can be enriched from mouse bone marrow using EasySep™ positive selection of c-kit<sup>+</sup> or Sca-1<sup>+</sup> cells.
  - No columns required.
  - High cell purity: utilizes the specificity of antibody-mediated selection.
- CFC are enriched over 20-fold with c-kit<sup>+</sup> selection and 10-fold with Sca-1<sup>+</sup> selection, with good recoveries (Table 2).
- EasySep™ selected c-kit<sup>+</sup> and Sca-1<sup>+</sup> cells are able to reconstitute multilineage hematopoiesis in irradiated recipients. The selected cells contain ~10 fold higher frequency of repopulating cells than the unseparated cells depending on the chosen protocol.

Together these observations demonstrate that EasySep™ positive selection of c-kit<sup>+</sup> and Sca-1<sup>+</sup> BM cells is a rapid and efficient means for the enrichment of primitive hematopoietic cells, including repopulating stem cells. This is a useful tool for further fractionation by FACS and for molecular and functional studies.