

## MoFlo™ Stem Cell Sorting

### Introduction

Purified populations of functional stem cells are of great interest to the biomedical community, both in the understanding of stem cell biology and in clinical transplant settings. Studies continue to define phenotypic markers, functional characteristics and *in vivo* reconstitutive activity for both hematopoietic and non-hematopoietic stem cells. Historically, the capture of these pluripotent cells has presented a significant laboratory challenge. In transplant programs, for instance, peripheral blood stem cells from a stimulated donor usually represent 0.01% – 1% of the collected white blood cells and bone marrow contains 0.5% – 5% stem cells.<sup>1</sup> A number of strategies have emerged for the purification of these cells, including high-throughput flow cytometry based on multiparametric immunophenotyping or immunophenotyping in combination with functional characteristics. The MoFlo High-Performance Cell Sorter has proved invaluable in these efforts. In the experiment shown here, the MoFlo was used to purify bone marrow side population cells.

### Materials and Methods

**Staining:** Bone marrow from tibias and femurs of 5- to 8-week-old C57BL/6 mice was flushed into HBSS in polypropylene centrifuge tubes. The nucleated cells were enumerated in this suspension, then pelleted and resuspended at  $1 \times 10^6$  cells/mL in pre-warmed DMEM. Hoechst 33342 was added to a final concentration of 5  $\mu$ g/mL.

Samples were mixed thoroughly and placed in a stable 37°C water bath for exactly 90 minutes, then immediately transferred to a 4°C centrifuge to pellet the cells. Cells were re-suspended in cold HBSS and maintained at 4°C. (If additional surface antibody labeling is desired, ensure that the cell suspension remains at 4°C. If desired, add propidium iodide at 2  $\mu$ g/mL to exclude non-viable cells from the sort.<sup>2</sup>)

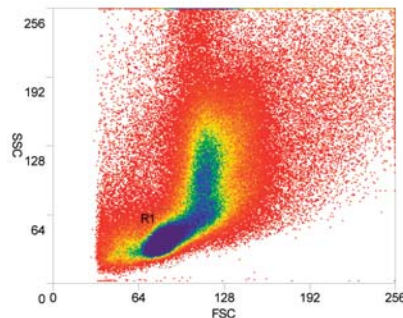
**Instrument Set-up.** A MoFlo with a 351 nm UV laser was configured to detect fluorescent emission both in the blue region, using a 450/20 bandpass filter, and in the red region, using a 675 eFLP filter.

### Results

Stained cells were placed on the MoFlo and a forward scatter vs. side scatter dot plot was used to gate the primary cell population (Figure 1a). A Hoechst Blue vs. Hoechst Red dot plot (Figure 1b) revealed the progenitor cells of interest, identified by their characteristic position to the left of the bulk cell population, i.e. the “side population”. These rare side population cells, a type of hematopoietic stem cell characterized by their ability to efflux Hoechst dye, were sorted for further investigation.

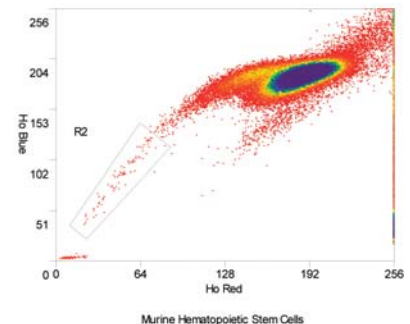
### Discussion

Purifying hematopoietic and non-hematopoietic stem cell populations,<sup>3-15</sup> including side populations,<sup>16-27</sup> has become routine with the MoFlo. The novel electronics, optics and fluidics of the MoFlo provide the power, precision and yield necessary to capture these rare events. Furthermore, the MoFlo uses a patented nozzle design to reduce turbulence and minimize the effects of acceleration on each cell. Thus, following purification, these cells are fully functional and capable of *in vivo* reconstitution, post-transplantation engraftment and long-term culture.<sup>28-34</sup>



**Figure 1a**

Plotting forward scatter vs. side scatter identifies the main cell population of interest.



**Figure 1b**

Following appropriate gating, the progenitor cells of interest are identified by their characteristic position to the left of the bulk cell population.

## References

- Ashcroft RG and Lopez PA. Commercial high-speed machines open new opportunities in high-throughput flow cytometry. *Journal of Immunological Methods* 2000; 243:13-24.
- Goodell MA. Hoechst 33342 HSC staining and stem cell purification protocol. Available at: [http://www.bcm.tmc.edu/genetherapy/goodell/new\\_site/index2.html](http://www.bcm.tmc.edu/genetherapy/goodell/new_site/index2.html)
- Akashi KX, He J, Chen H, Iwasaki C, Niu B, Steenhard J et al. Transcriptional accessibility for genes of multiple tissues and hematopoietic lineages is hierarchically controlled during early hematopoiesis. *Blood* 2003; 101:383.
- Goan SR, Junghahn I, Wissler M, Becker M, Aumann J, Just U et al. Donor stromal cells from human blood engraft in NOD/SCID mice. *Blood* 2002; 96:3971.
- Habibian HK, Peters SO, Hsieh CC, Wu J, Vergilis K, Grimaldi CI et al. The fluctuating phenotype of the lymphohematopoietic stem cell with cell cycle transit. *J Exp Med* 1998; 188:393.
- Henckaerts E, Geiger H, Langer JC, Rebollo P, Van Zant G, Snoeck HW. Genetically determined variation in the number of phenotypically defined hematopoietic progenitor and stem cells and in their response to early-acting cytokines. *Blood* 2002; 99:3947.
- Jackson KA, Mi T, Goodell MA. Hematopoietic potential of stem cells isolated from murine skeletal muscle. *PNAS* 1999; 96:14482.
- Kirby S, Walton W, Smithies O. Hematopoietic stem cells with controllable tEpoR transgenes have a competitive advantage in bone marrow transplantation. *Blood* 2000; 95: 3710.
- Kuehnle I, Huls MH, Liu Z, Semmelmann M, Krance RA, Brenner MK et al. CD20 monoclonal antibody (rituximab) for therapy of Epstein-Barr virus lymphoma after hematopoietic stem-cell transplantation. *Blood* 2000; 95:1502.
- Manz MG, Miyamoto T, Akashi K, Weissman IL. Prospective isolation of human clonogenic common myeloid progenitors. *PNAS* 2002; 99:11872.
- McKinney-Freeman SL, Jackson KA, Camargo FD, Ferrari G, Mavilio F, Goodell MA. Muscle-derived hematopoietic stem cells are hematopoietic in origin. *PNAS* 2002; 99:1341.
- Okuno Y, Huettner CS, Radomska HS, Petkova V, Iwasaki H, Akashi K et al. Distal elements are critical for human CD34 expression in vivo. *Blood* 2002; 100:4420.
- Schaniel C, Bruno L, Melchers F, Rolink AG. Multiple hematopoietic cell lineages develop in vivo from transplanted Pax5-deficient pre-B1-cell clones. *Blood* 2002; 99:472.
- Wright DE, Bowman EP, Wagers AJ, Butcher EC, Weissman IL. Hematopoietic stem cells are uniquely selective in their migratory response to chemokines. *J Exp Med* 2002; 195:1145.
- Zhong JF, Zhan Y, Anderson WF, Zhao Y. Murine hematopoietic stem cell distribution and proliferation in ablated and nonablated bone marrow transplantation. *Blood* 2002; 100:3521.
- Asakura A, Seale P, Girgis-Gabardo A, Rudnicki MA. Myogenic specification of side population cells in skeletal muscle. *J Cell Biol* 2002; 159:123-134.
- Falciatori I, Borsellino G, Haliassos N, Boitani C, Corallini S, Battistini L et al. Identification and enrichment of spermatogonial stem cells displaying side-population phenotype in immature mouse testis. *FASEB J* 2003; The FASEB Journal Express Article doi:10.1096/fj.03-0744fje; Published online December 19, 2003.
- Jackson KA, Majka SM, Wang H, Pocius J, Hartley CJ, Majesky MW, et al. Regeneration of ischemic cardiac muscle and vascular endothelium by adult stem cells. *J Clin Invest* 2001; 107: 1395-402.
- Klarmann K, Ortiz M, Davies M, Keller JR. identification of in vitro growth conditions for c-Kit-negative hematopoietic stem cells. *Blood* 2003; 102: 3120-3128.
- Oh H, Bradfute SB, Gallardo TD, Nakamura T, Gausin V, Mishina Y et al. Cardiac progenitor cells from adult myocardium: homing, differentiation, and fusion after infarction. *PNAS* 2003; 21: 12313-12318.
- Okuno Y, Iwasaki H, Huettner CS, Radomska HS, Gonzalez DA, Tenen DG et al. Differential regulation of the human and murine CD34 genes in hematopoietic stem cells. *PNAS* 2002; 99: 6246-51.
- Majka SM, Jackson KA, Kienstra KA, Majesky MW, Goodell MA, Hirschi KK. Distinct progenitor populations in skeletal muscle are bone marrow derived and exhibit different cell fates during vascular regeneration. *J Clin Invest* 2003; 111: 71-79.
- Mogi M, Yang J, Lambert JF, Colvin GA, Shiojima I, Skurk C et al. Akt Signaling Regulates Side Population Cell Phenotype via Bcrp1 Translocation. *J Biol Chem* 2003; 278: 39068-39075.
- Pearce DJ, Ridler CM, Simpson C, Bonnet D. Multi-parameter analysis of murine bone marrow side population cells. *Blood* 2003; *Blood First Edition Paper*, prepublished online Nov 26 2003, DOI 10.1182/Blood-2003-09-3281.
- Radomska HS, Gonzalez DS, Okuno Y, Iwasaki H, Nagy A, Akashi K et al. Transgenic targeting with regulatory elements of the human CD34 gene. *Blood* 2002; 100: 4410-4419.
- Summer R, Kotton DN, Sun X, Ma B, Fitzsimmons K, Fine A. Side population cells and Bcrp1 expression in lung. *American Journal of Physiology, Lung Cellular and Molecular Physiology* 2003; 285: 97-104.
- Wulf GG, Wang RY, Kuehnle I, Weidner D, Marini F, Brenner MK et al. A leukemic stem cell with intrinsic drug efflux capacity in acute myeloid leukemia. *Blood* 2001; 98: 1166-73.
- Chatterjee S et al. Transduction of primitive human marrow and cord blood-derived hematopoietic progenitor cells with adeno-associated virus vectors. *Blood* 1999; 93(6): 1882-1894.
- Hulspas R et al. Characterization of neurosphere cell phenotypes by flow cytometry. *Cytometry* 2000; 40: 245-250.
- Shih C et al. Long-term ex vivo maintenance and expansion of transplantable human hematopoietic stem cells. *Blood* 1999; 94(5): 1623-1636.
- Shih CC, Hu MC, Hu J, Weng Y, Yazaki PJ, Medeiros J et al. A secreted and LIF-mediated stromal cell-derived activity that promotes ex vivo expansion of human hematopoietic stem cells. *Blood* 2000; 95:1957.
- Shih CC, Weng Y, Mamelak A, LeBon T, Hu MC, Forman SJ. Identification of a candidate human neurohematopoietic stem-cell population. *Blood* 2001; 98:2412.
- Spyridonidis A et al. Purging of mammary carcinoma cells during ex vivo culture of CD34+ hematopoietic progenitor cells with recombinant immunotoxins. *Blood* 1998; 91(5): 1820-1827.
- Stewart FM et al. Lymphohematopoietic engraftment in minimally myeloablated hosts. *Blood* 1998; 91(10): 3681-3687.

## Acknowledgements

Special thanks to Karen Helm (University of Colorado Health Sciences Center) and Margaret Goodell, PhD (Center for Cell & Gene Therapy, Baylor College of Medicine) for data and protocols used in this report.

For Research Use Only. Not for use in Diagnostic Procedures.

Australia, Gladesville (61) 2 9844-6000 Canada, Mississauga (1) 905 819 1234 China, Beijing (86) 10 6515 6028  
Czech Republic, Prague (420) 272 01 73 32 Eastern Europe, Middle East, North Africa, South West Asia: Switzerland, Nyon (41) 22 365 3707  
France, Villepinte (33) 1 49 90 90 00 Germany, Krefeld (49) 2151 33 35 Hong Kong (852) 2814 7431  
India, Mumbai (91) 22 3080 5101 Italy, Cassina de' Pecchi, Milan (39) 02 953921 Japan, Tokyo (81) 3 5530 8500 Korea, Seoul (82) 2 404 2146  
Latin America (1) (305) 380 4709 Mexico, Mexico City (001) 52 55 9183 2800 Netherlands, Mijdrecht (31) 297 230630 Puerto Rico (787) 747 3335  
Singapore (65) 6339 3633 South Africa/Sub-Saharan Africa, Johannesburg (27) 11 564 3203 Spain, Madrid (34) 91 3836080  
Sweden, Bromma (46) 8 564 85 900 Switzerland, Nyon (41) 0800 850 810 Taiwan, Taipei (886) 2 2378 3456 Turkey, Istanbul (90) 216 309 1900  
UK, High Wycombe (44) 01494 441181 USA, Fullerton, CA (1) 800 742 2345

