

EasySep™ Kits for Immunomagnetic Purification of CD138+ cells from Mouse and Rat Samples

Trevor Rogers¹, Kathleen MacDonald¹, Whalley Fong¹, Peter Repenning¹, Alexisann Maxwell¹, Mark Brown¹, Terry Thomas¹, and Allen Eaves^{1,2}

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

trevor.rogers@stemcell.com

Introduction

CD138 is found on several B cell subsets but is highly expressed on antibody-secreting cells (plasma cells and plasmablasts). Because CD138+ antibody-secreting cells represent a small fraction of the total B cell population (0.4 ± 0.1% (n = 52), 1.6 ± 1.1% (n = 43), and 0.6 ± 0.5% (n = 18) when using naïve C57BL/6 mice, immunized BALB/c mice, or naïve rat (Sprague Dawley and Wistar) spleens, respectively), pre-enrichment of CD138+ cells can improve downstream assays and processes such as hybridoma generation and screening. Therefore, we have developed a facile immunomagnetic method to isolate CD138+ cells from mouse and rat samples.

The column-free immunomagnetic method allows for isolation of CD138+ cells in less than 25 minutes from mouse or rat spleen, bone marrow, and other tissues using fast and simple EasySep™ technology. Bispecific antibody complexes are used to cross-link CD138+ cells to dextran-coated magnetic particles and the labeled cells are separated from the unwanted cells using an EasySep™ magnet. Single-cell suspensions of splenocytes and bone marrow cells do not require red blood cell lysis prior to isolation.

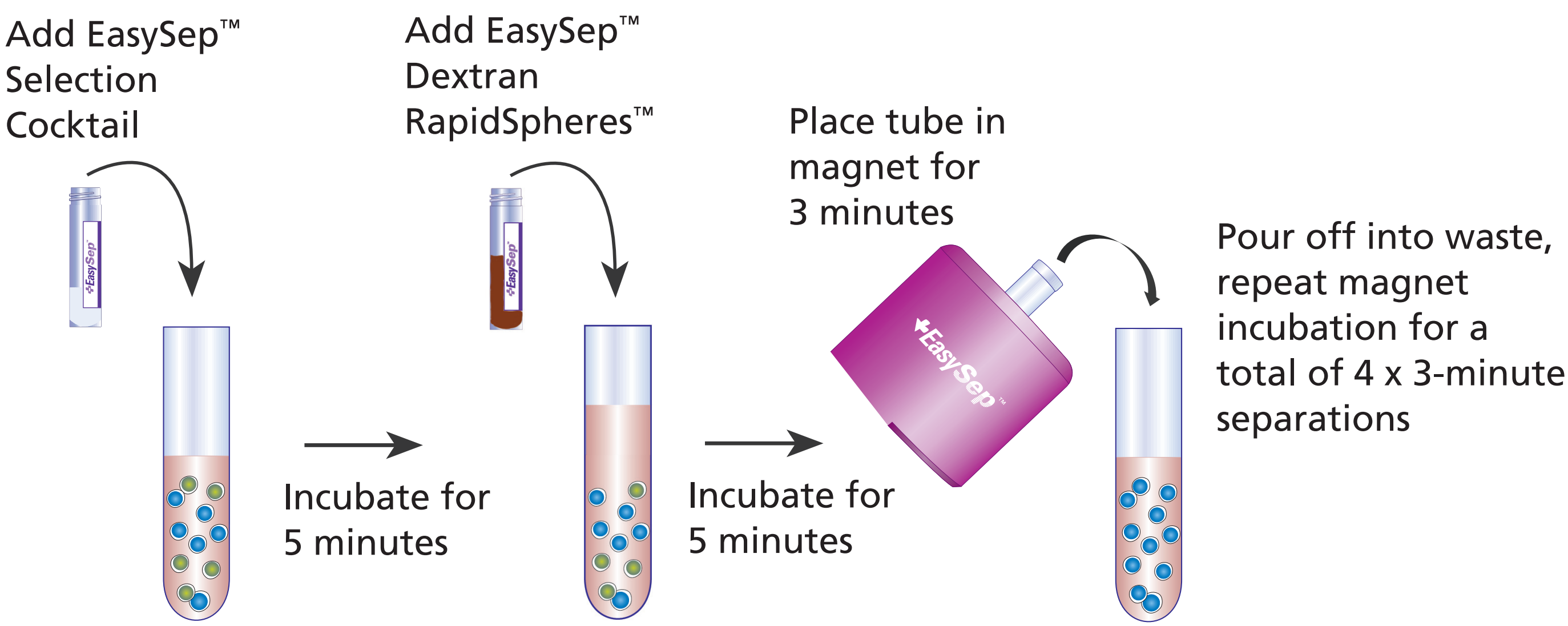
Using this method, the total viable CD138+ content of the isolated fraction is typically 75.0 ± 10.2% (n = 41), 77.9 ± 1.6% (n = 4), and 55.5 ± 10.3% (n = 52), when using naïve C57BL/6 mice, immunized BALB/c mice, or naïve rat spleens, respectively, while the high-expressing CD138+ antibody-secreting cell content of the isolated fraction is typically 49.5 ± 16.2% (n = 41), 71.6 ± 3.4% (n = 4), and 48.6 ± 10.8% (n = 52), respectively. The isolated CD138+ populations remain functional and are immediately ready for downstream applications, including antibody development projects.

Methods

Preparation of Starting Cell Suspension

To prepare a single-cell suspension of splenocytes, spleens from naïve C57BL/6 mice, immunized BALB/c mice, or naïve Sprague Dawley or Wistar rats were disrupted in phosphate-buffered saline (PBS) containing 2% fetal bovine serum (FBS) and 1 mM EDTA. To remove aggregates and debris, cells were passed through a 70 µm mesh nylon strainer and centrifuged at 120 x g for 10 minutes. A mononuclear cell suspension was prepared from rat bone marrow by flushing cells from the femur and tibia bones into PBS/2% FBS/1 mM EDTA using a 23-gauge needle. All cells were resuspended at 1 x 10⁸ cells/mL in PBS/2% FBS/1 mM EDTA prior to EasySep™ separation.

Figure 1. EasySep™ protocol for isolating mouse or rat CD138+ cells from splenocytes or bone marrow suspensions



The procedure involves the addition of the EasySep™ Cocktail and EasySep™ Dextran RapidSpheres™ to the cells, followed by immunomagnetic separation.

Results

Table 1. Purities and average yields of CD138+ cells isolated from rat spleen or bone marrow cells by EasySep™

CD138 Total					
Sample	Start		Isolated		
	n	Purity (% ± SD)	n	Purity (% ± SD)	Target cells recovered per 1 x 10 ⁸ start cells
Naïve Rat Spleen	17	0.9 ± 0.8	52	55.5 ± 10.3	0.39 x 10 ⁶
Naïve Rat Bone Marrow	15	8.8 ± 5.1	87	62.7 ± 9.0	1.30 x 10 ⁶
Plasma cells/Plasmablasts					
Sample	Start		Isolated		
	n	Purity (% ± SD)	n	Purity (% ± SD)	Target cells recovered per 1 x 10 ⁸ start cells
Naïve Rat Spleen	18	0.6 ± 0.5	52	48.6 ± 10.8	0.28 x 10 ⁶
Naïve Rat Bone Marrow	15	0.4 ± 0.3	86	9.2 ± 7.8	0.10 x 10 ⁶

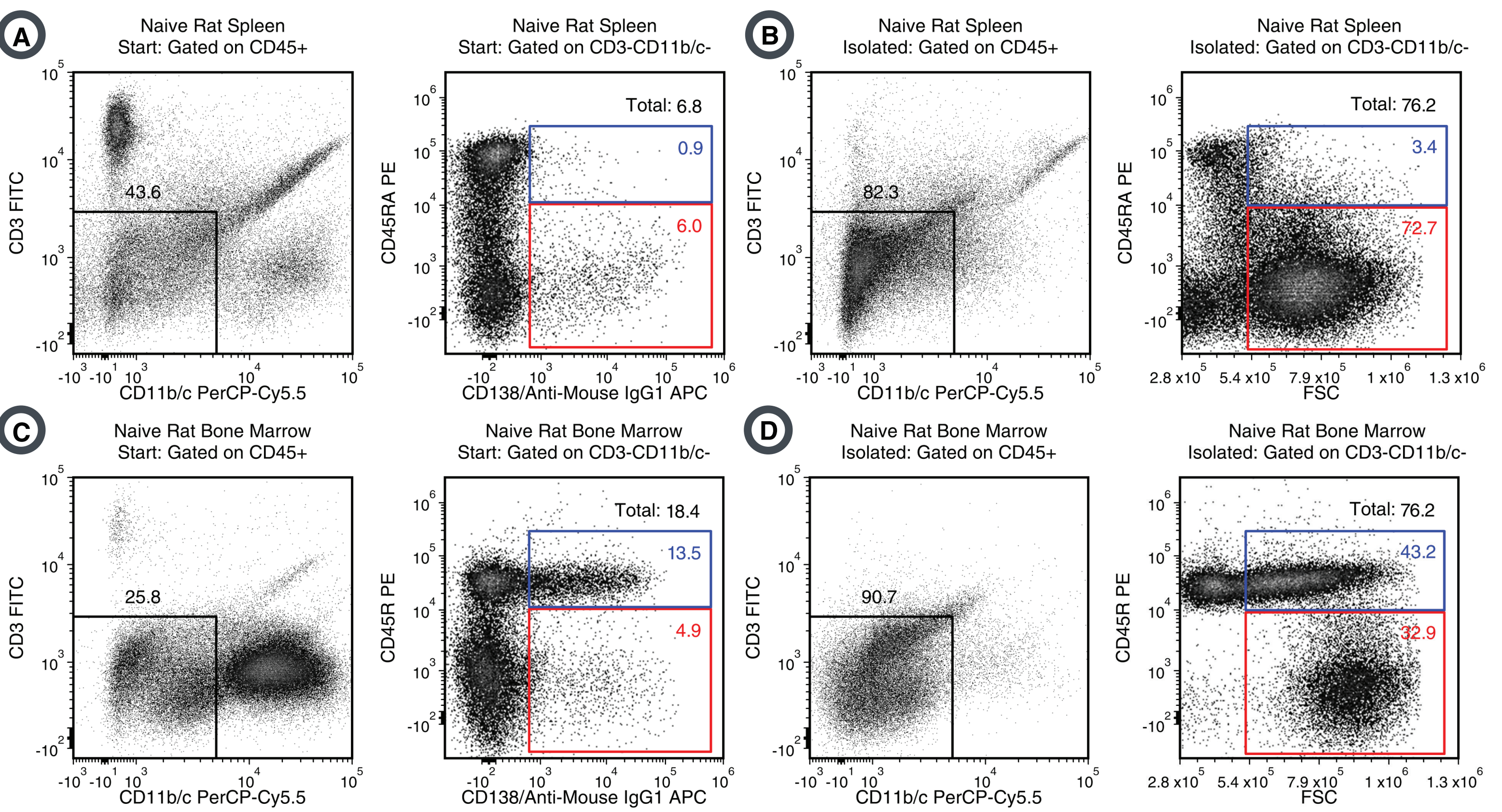
Purities and average cell yields of isolated rat CD138+ cells are based on the percentage of viable CD45+CD3+CD11b/c- cells. Purity values are expressed as mean ± SD. Average cell yield is presented as the number of cells recovered per 1 x 10⁸ start cells. Results include data for EasySep™ Magnet, "The Big Easy" Magnet, and EasyEights™ Magnet.

Table 2. Purities and average cell yields of mouse CD138+ cells isolated by EasySep™

CD138 Total					
Sample	Start		Isolated		
	n	Purity (% ± SD)	n	Purity (% ± SD)	Target cells recovered per 1 x 10 ⁸ start cells
Naïve Mouse Spleen	52	6.0 ± 2.1	41	75.0 ± 10.2	0.36 x 10 ⁶
Immunized Mouse Spleen	43	5.9 ± 2.3	4	77.9 ± 1.6	2.48 x 10 ⁶
Plasma cells/Plasmablasts					
Sample	Start		Isolated		
	n	Purity (% ± SD)	n	Purity (% ± SD)	Target cells recovered per 1 x 10 ⁸ start cells
Naïve Mouse Spleen	52	0.4 ± 0.1	41	49.5 ± 16.2	0.24 x 10 ⁶
Immunized Mouse Spleen	43	1.6 ± 1.1	4	71.6 ± 3.4	1.47 x 10 ⁶

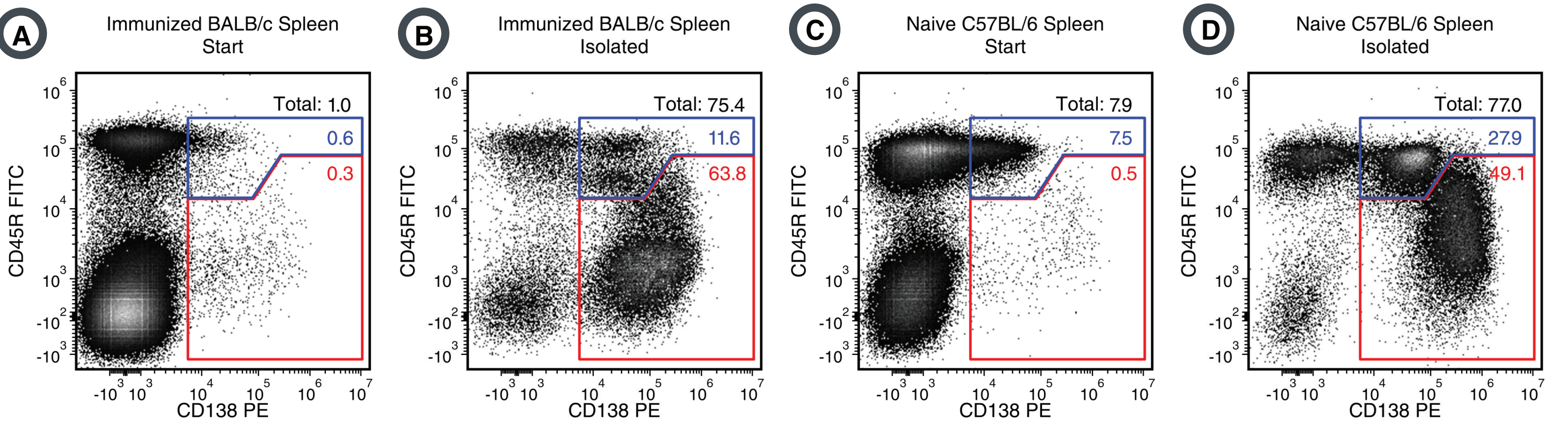
Purities and average cell yields of isolated mouse CD138+ cells is based on percentage of viable cells. Average cell yield is indicated as the number of cells recovered per 1 x 10⁸ start cells from naïve C57BL/6 or immunized BALB/c mice spleens (purity values are expressed as mean ± SD). Results include data for EasySep™ Magnet, "The Big Easy" Magnet, and EasyEights™ Magnet.

Figure 2. Representative flow cytometry data before and after EasySep™ isolation of CD138+ cells from naïve rat spleen or bone marrow samples



The viable CD138+ content (gated on CD45+CD3+CD11b/c- cells) is shown for start (A) and isolated (B) splenocytes, and for start (C) and isolated (D) bone marrow cells from naïve Sprague Dawley rats. Cells were assessed by flow cytometry with antibodies against CD45, CD3, CD11b/c, and either CD45RA (splenocytes) or CD45R (bone marrow cells) in the presence of DAPI. Start CD138+ cells were identified with an anti-rat CD138 antibody, followed by a rat anti-mouse IgG1 antibody, APC. Isolated plasma cells and blasts are identified as large FSC/CD45RA^{Low/-} (spleen) or large FSC/CD45R^{Low/-} (bone marrow).

Figure 3. Representative flow cytometry data before and after EasySep™ isolation of CD138+ cells from the spleens of immunized or naïve mice



The viable CD138+ content is shown for start (A) and isolated (B) cells from immunized BALB/c mice, and for start (C) and isolated (D) cells from naïve C57BL/6 mice. Samples were assessed by flow cytometry using antibodies against CD138 and CD45R in the presence of DAPI. Plasma cells/plasmablasts are identified as CD138^{High} and CD45R^{Low/-}.

Summary

- Fast, easy-to-use, and column-free isolation of CD138+ cells directly from mouse or rat tissues in as little as 25 minutes
- No red blood cell lysis required prior to isolation
- Isolated CD138+ populations remain functional and are immediately ready for downstream assays and processes such as hybridoma generation and screening