

# A Rapid Method to Enrich Specific Lymphocyte Populations (T cells, B cells, Total Lymphocytes) from Whole Blood

Karina L. McQueen<sup>1</sup>, Jenna L. Warren<sup>1</sup>, Allen C. Eaves<sup>1,2</sup> and Terry E. Thomas<sup>1</sup>

<sup>1</sup>StemCell Technologies Inc., 570 West 7th Ave., Vancouver, BC, V5Z 1B3, Canada <sup>2</sup>Terry Fox Laboratory, 675 West 10th Ave, Vancouver, BC, V5Z 1L3, Canada

## Summary

The isolation of specific lymphocyte populations is essential for many HLA applications. While density centrifugation with Ficoll™ is used by most laboratories, it is a lengthy process and can result in lymphocyte loss of up to 50%. Isolation of specific cell types directly from whole blood requires a red blood cell lysis/wash step which is time consuming and is often toxic to certain cell subsets.

We have developed a rapid (20 minute) immunomagnetic cell separation system (EasySep®) that enriches for T cells, B cells or total lymphocytes from whole blood. This method does not require Ficoll™ density centrifugation or lysis to remove the red blood cell burden from the sample, but instead relies on hetastarch sedimentation (HetaSep™) which is less toxic, produces a high yield of lymphocytes and typically can be performed in 15 minutes.

Assessment by flow cytometry yields average purities greater than 90% for all cell types. As the cells of interest are not labeled with antibody, they are immediately available for all downstream HLA applications.

## Methods

### Preparation of sample using HetaSep™:

Whole blood was collected in a blood collection tube containing heparin or ACD. One part HetaSep™ (Catalog #07806) was added to 5 parts whole blood and mixed. The sample was placed in a 37°C incubator and allowed to settle until the red blood cell interface was at approximately 40% of the total volume. The supernatant was then harvested, washed once and centrifuged at room temperature at 120 x g for 10 minutes with the brake off. Cells were resuspended in 1/10th the original starting volume of whole blood.

### The starting cell number:

From a starting blood volume of 5.0 mL, the number of nucleated cells used per experiment ranged from 7.5x10<sup>6</sup> – 3.0x10<sup>7</sup>.

### EasySep® enrichment kits used to isolate specific lymphocyte subsets from whole blood:

Total Lymphocyte Enrichment: Catalog #19961HLA

T Cell Enrichment: Catalog # 19951HLA

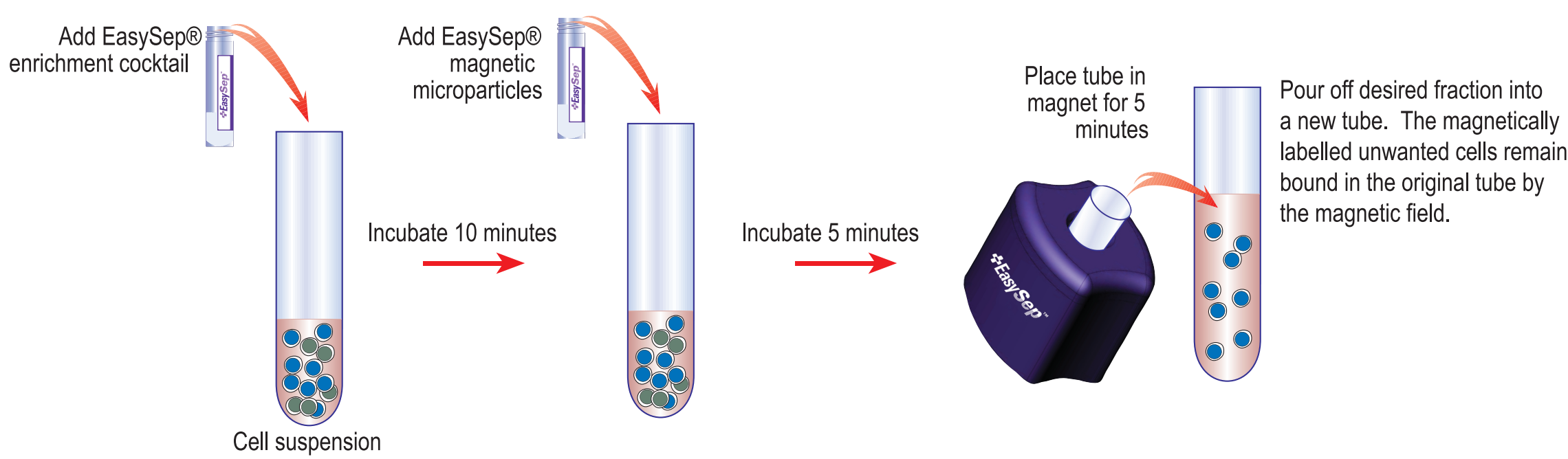
B Cell Enrichment: Catalog # 19954HLA

Following HetaSep™ treatment of the whole blood, cells are labeled with a cocktail of antibodies targeting unwanted cells. Cells are then coupled to magnetic particles and the sample is placed in a magnet. Labeled unwanted cells remain in the sample tube in the magnet, while the unlabeled cells of interest are removed to a new tube.

Cell isolation can be performed manually, or can be automated using RoboSep®.

Purity of selected cell populations was determined by flow cytometry. Total lymphocytes were defined as CD3<sup>+</sup>/CD19<sup>+</sup>; T cells were defined as CD3<sup>+</sup>; B cells were defined as CD19<sup>+</sup>. Only nucleated cells (CD45<sup>+</sup>) were included in this assessment.

Figure 1: Schematic Representation of the Manual EasySep® Procedure.



## Results

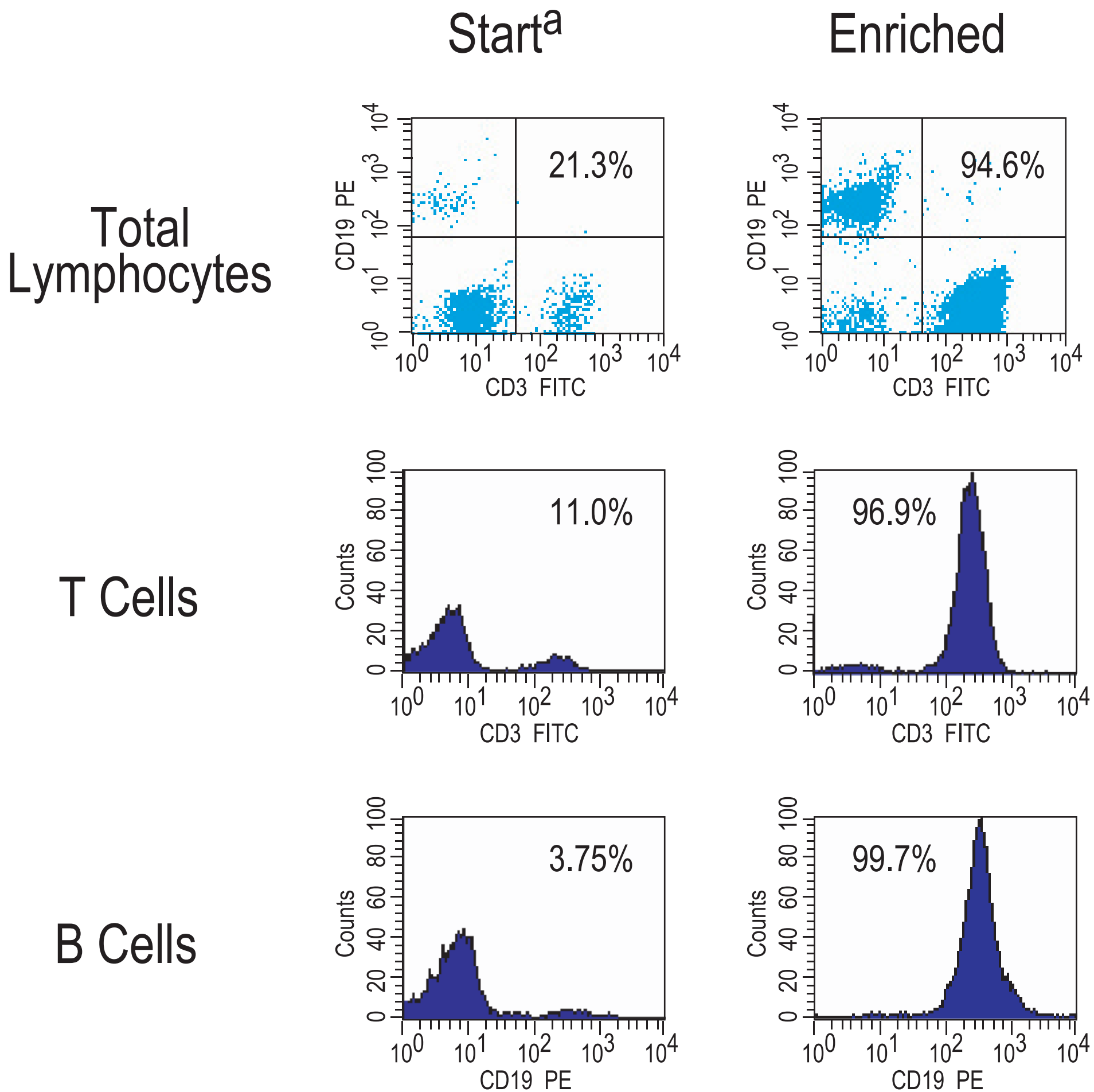
Table 1. Percent purity, recovery and average number of enriched cells obtained from 5.0 mL of whole blood using the EasySep® Total Lymphocyte (TL), T cell and B cell enrichment kits. Purities were determined by flow cytometry. All samples were gated on CD45<sup>+</sup>, viable (PI negative) cells. Values are expressed as means +/- 1 standard deviation.

	n	% Start <sup>a</sup>	% Purity	% Recovery <sup>a</sup>	# Enriched Cells (x10 <sup>6</sup> ) <sup>b</sup>
TL	8	26.9 ± 8.8	93.6 ± 2.3	73.4 ± 12.6	2.6
T cell	16	21.6 ± 8.6	96.6 ± 1.6	76.8 ± 25.5	2.2
B cell	17	4.1 ± 1.6	96 ± 5.6	58.5 ± 20.3	0.4

<sup>a</sup> From HetaSep™ - treated whole blood

<sup>b</sup> From 5.0 mL of whole blood starting sample

Figure 2. Typical FACS plots before and after enrichment of selected cells (gated on CD45<sup>+</sup> cells).



<sup>a</sup> From HetaSep™-treated whole blood

Table 2. Direct comparison between the number of purified cells obtained from 10 mL of whole blood using either EasySep® enrichment or a competitor's negative selection system. EasySep® HLA enrichment was performed on HetaSep™-treated blood, whereas the competitor's enrichment was performed on a mononuclear cell suspension. In both systems, 10 mL of whole blood from the same donor was used as the starting sample.

Cell Type	Average Number of Enriched Cells (x10 <sup>6</sup> )		
	EasySep®	Competitor	Fold Difference
T Cells <sup>a</sup>	6.5	3.5	1.8
B Cells <sup>b</sup>	0.9	0.43	2.6

<sup>a</sup> average from 3 experiments

<sup>b</sup> average from 6 experiments

## Conclusions

- Total lymphocytes, T cells or B cells can be rapidly isolated in approximately 20 minutes
- The red blood cell burden of the sample is reduced prior to cell isolation by hetastarch sedimentation using HetaSep™
- No layering of Ficoll™ or post-enrichment lysis of the sample is required
- Purity of total lymphocytes ranges from 90.2 – 96.9%, purity of T cells ranges from 93.1 – 98%, while purity of B cells ranges from 81.5 – 99.4%
- A 5.0 mL whole blood start sample yields enough cells for many HLA applications
- Purified cells are unlabeled (no antibodies or magnetic particles) and are immediately available for all HLA applications



StemCell Technologies Inc

The Cell Experts™ | www.stemcell.com

In North America  
Tel: 1.604.877.0713  
Fax: 1.604.877.0704  
Toll Free Tel: 1.800.667.0322  
Toll Free Fax: 1.800.567.2899  
e-mail: info@stemcell.com

In Europe  
Tel: +33.(0).4.76.04.75.30  
Fax: +33.(0).4.76.18.99.63  
e-mail: info@stemcellfrance.com

In the United Kingdom  
Tel: +44.(0).20.7537.7565  
Fax: +44.(0).20.7515.5408  
Toll Free within UK:  
Tel: 0800.731.27.14  
Fax: 0800.731.27.13  
e-mail: info@stemcellgb.com