

# A simple and rapid method for the enrichment of mouse naïve CD4<sup>+</sup> T cells from spleen

Vesna Posarac, Armon Molavi, Karina McQueen, and Terry E. Thomas  
STEMCELL Technologies Inc., Vancouver, BC, Canada

## Abstract

Naïve CD4<sup>+</sup> T cells are a mature subset of CD4<sup>+</sup> T cells with no previous antigen exposure. The CD62L<sup>high</sup> CD44<sup>low</sup> naïve phenotype cells circulate throughout the secondary lymphoid organs where they become activated by foreign antigens presented on MHC class II molecules. Activation is marked by phenotypic changes (down- and up-regulation of CD62L and CD44, respectively), proliferation, and acquisition of effector functions. Current protocols for the isolation of naïve CD4<sup>+</sup> T cells are time-consuming and require the use of columns. We have developed a one-step, column-free, immunomagnetic cell separation (EasySep™) method for the isolation of naïve CD4<sup>+</sup> T cells from single-cell suspensions of splenocytes. Non-CD4<sup>+</sup> T cells, T regulatory cells, and memory CD4<sup>+</sup> T cells are targeted for depletion using biotinylated antibodies cross-linked to streptavidin-coated magnetic particles. The labeled cells are separated using an EasySep™ magnet and the desired naïve CD4<sup>+</sup> T cell fraction is poured off. The entire protocol is performed in 15 minutes and can be fully automated using RoboSep™. The average purities and recoveries of CD4<sup>+</sup>CD62L<sup>high</sup> CD44<sup>low</sup> cells are 93.1 ± 2.1% and 28.8 ± 7.6%, respectively (n=22). This new method will be invaluable to researchers studying T cell differentiation, signaling pathways, and immune response to infectious disease.

## Methods

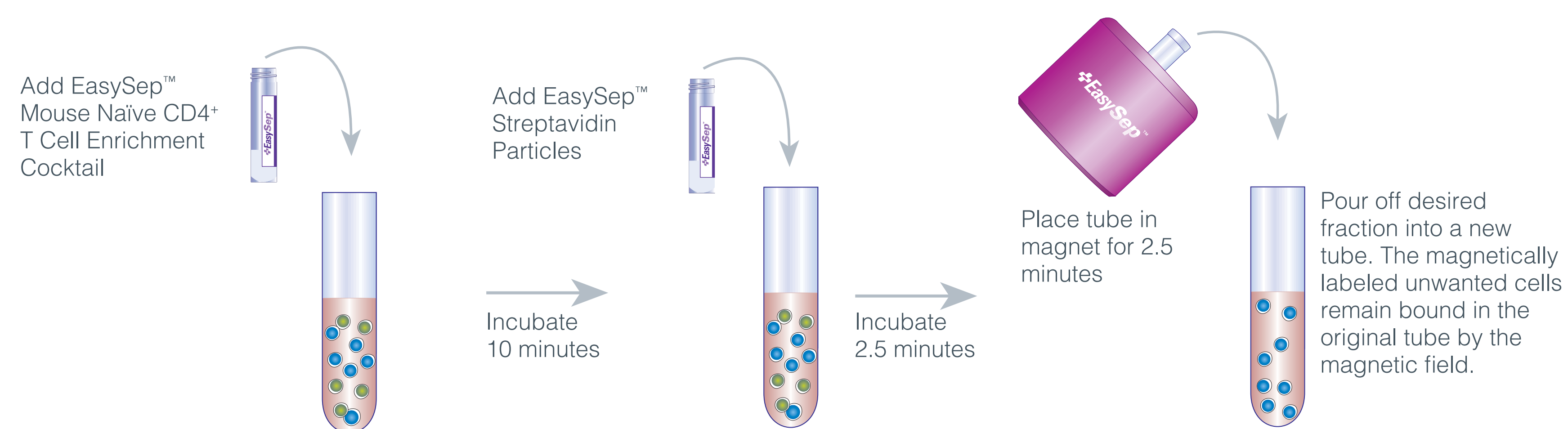
### Preparation of Starting Cell Suspension

To prepare a single-cell suspension, spleens were disrupted in phosphate buffered saline (PBS) + 2% fetal bovine serum (FBS). The cells were centrifuged at 300 x g for 10 minutes and resuspended at 1x10<sup>8</sup> cells per ml in PBS + 2% FBS with 5% normal rat serum.

### EasySep™ Labeling of Mouse Cells

Unwanted cells, including CD4<sup>-</sup> cells, T regulatory cells, and CD44<sup>high</sup> CD4<sup>+</sup> T cells are specifically labeled with biotinylated antibodies cross-linked to streptavidin coated magnetic particles. The magnetically labeled cells are then separated from unlabeled cells using an EasySep™ magnet (Figure 1), and the desired naïve CD4<sup>+</sup> T cell fraction is poured off.

**FIGURE 1: EasySep™ procedure for column-free enrichment of naïve CD4<sup>+</sup> T cells from mouse splenocytes**



This procedure can be fully automated using RoboSep™.

### Purity Assessment

EasySep™-isolated naïve CD4<sup>+</sup> T cells were assessed by flow cytometry after staining with CD44 FITC, CD62L PE, CD4 APC, and 7-AAD. Naïve CD4<sup>+</sup> T cells are CD4<sup>+</sup>CD62L<sup>high</sup>CD44<sup>low</sup> (Figure 2).

## Results

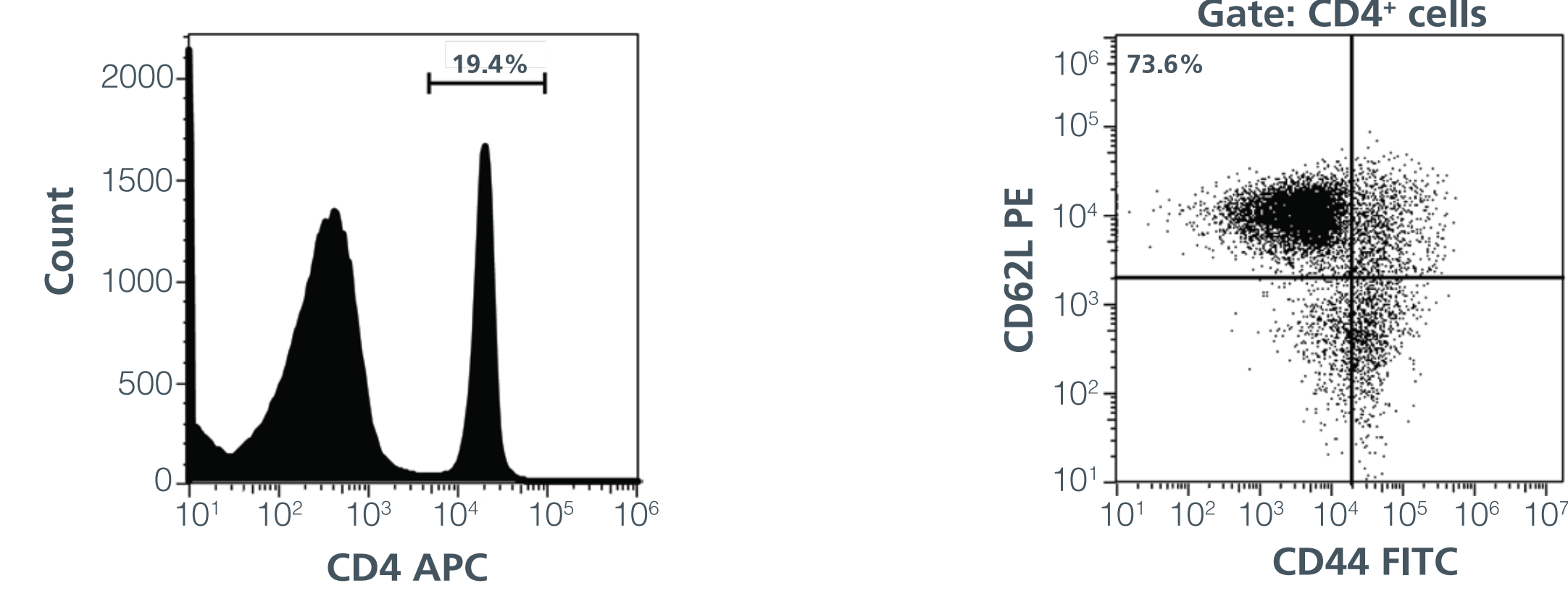
**TABLE 1: Purity and recovery of naïve CD4<sup>+</sup> T cells enriched from mouse splenocytes by EasySep™ or RoboSep™**

Method	n	% Purity	% Recovery
EasySep™	16	93.8 ± 1.6	29.6 ± 7.8
RoboSep™	6	91.2 ± 2.4	26.8 ± 7.5

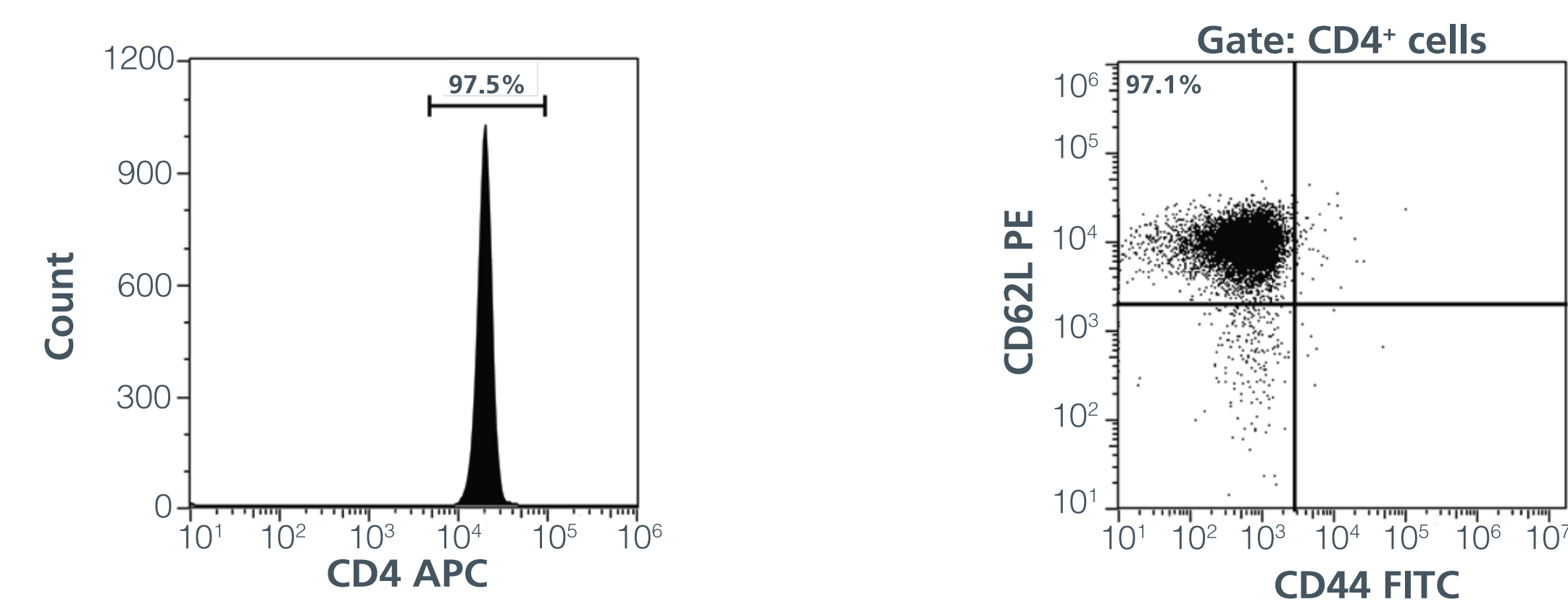
Purities were determined by flow cytometry. All samples gated on viable (7-AAD<sup>-</sup>) cells. Naïve CD4<sup>+</sup> T cells are defined as CD4<sup>+</sup>CD62L<sup>high</sup>CD44<sup>low</sup>. Values are expressed as means ± 1 SD.

**FIGURE 2: Flow cytometric assessment of naïve CD4<sup>+</sup> T cells before and after enrichment using EasySep™**

**Start:** 14.3% CD4<sup>+</sup>CD62L<sup>high</sup>CD44<sup>low</sup> viable cells



**Enriched:** 94.7% CD4<sup>+</sup>CD62L<sup>high</sup>CD44<sup>low</sup> viable cells



**TABLE 2: Comparison of naïve CD4<sup>+</sup> T cell isolation protocols using EasySep™/RoboSep™ or the column-based competitor kit**

	EasySep™	RoboSep™	Competitor
Total time	15 min	22 min	1 hr 50 min
Columns	0	0	2
Centrifugations	0	0	4
Isolation method	untouched	untouched	positive selection

## Conclusions

- Isolate naïve CD4<sup>+</sup> T cells from mouse splenocytes in 15 minutes.
- Naïve CD4<sup>+</sup> T cell isolation can be fully automated with RoboSep™.
- Average purities and recoveries for naïve CD4<sup>+</sup> T cell enrichments are 93.1% ± 2.1% and 28.8% ± 7.6%, respectively.