

Quick and Easy Isolation of Immune Cells From Large-Volume Samples

Vesna Posarac¹, G Neil MacDonald¹, Chris A. Buck¹, Eric Toombs¹, Savannah D. Gellner¹, Hayley Reeves¹, Susan de Jong¹, Mark E. Williamson¹, Oliver Egeler¹, Bob Dalton¹, Andy I. Kokaji¹, Allen C. Eaves^{1,2}, Sharon A. Louis¹, and Frann Antignano¹

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

ABSTRACT

Large-scale cell isolation is commonly performed in labs and core facilities for cell banking and drug discovery, and as a critical step in cell therapy manufacturing. Current methods can be a significant bottleneck in a lab's workflow, often requiring a full day for sample processing and cell isolation. To address this need, we have developed two methods: manual column-free isolation using the Easy 250 EasySep™ magnet, and automated cell isolation in a closed system using RoboSep™-C.

The Easy 250 magnet allows users to isolate cells from a full leukopak in under 30 minutes by simply pipetting out their target cells. Negative selection protocols have been optimized to achieve 95.3% T cell, 97.2% CD4⁺ T cell, 91.9% CD8⁺ T cell, 99.4% B cell, 96.5% NK cell, and 91.5% monocyte purities. Positive selection protocols obtain 95.4% CD3⁺ cell, 94.4% CD4⁺ T cell, 93.9% CD8⁺ T cell, and 96.2% CD14⁺ cell purities.

RoboSep™-C automates this cell isolation procedure, along with the cell washing steps for sample preparation, in as little as 50 minutes. The instrument features a scale tower, pump and clamp modules to direct fluid through defined paths, and a magnet for separation of labeled cells. The system uses a sterile single-use tubing set that incorporates a cartridge for cell washing and concentration, and a magnet chamber for cell separation. Starting with fresh leukopaks, we obtained 95.9% T cell, 95.8% CD4⁺ T cell, 89.6% CD8⁺ T cell, 92.6% monocyte, and 90.5% NK cell purities following negative selection.

These approaches offer efficient and user-friendly cell isolation that allows researchers to scale up their operations, and can be easily integrated upstream of existing workflows.

METHODS

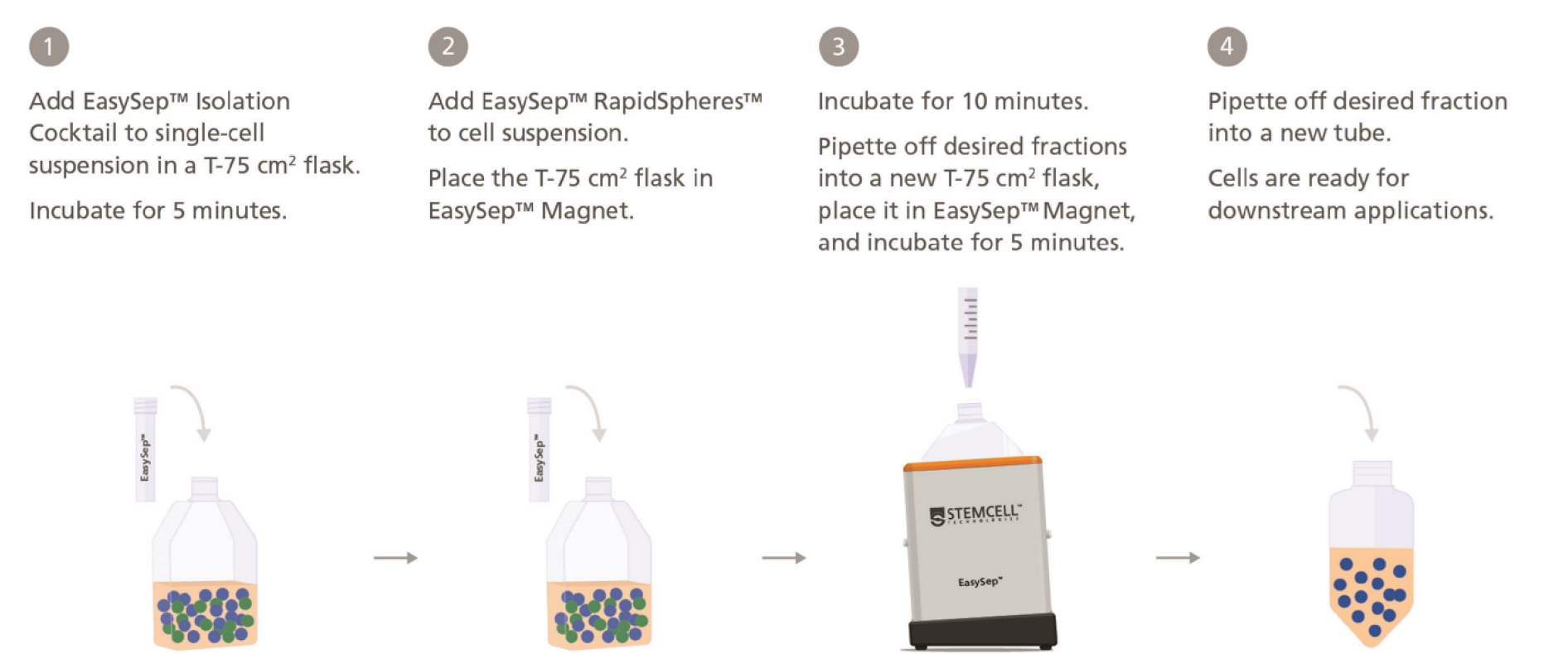


FIGURE 1. Typical Cell Isolation Protocol Using the Easy 250 EasySep™ Magnet.

To prepare starting samples for cell isolation, leukopaks are washed by diluting with an equal volume of EasySep™ Buffer and centrifuged at 300 x *g* for 10 minutes. Optionally, red blood cell (RBC) lysis can be performed using ammonium chloride. Washed cells are resuspended at 5 x 10⁷ cells/mL and up to 225 mL for negative selection, or 1 x 10⁸ cells/mL and up to 125 mL for positive selection. Unwanted cells (negative selection) or target cells (positive selection) are labeled using antibody complexes and magnetic particles and are separated in a T-75 cm² flask using the Easy 250 EasySep™ Magnet. Estimated time for the cell isolation protocol is 20 - 30 minutes.

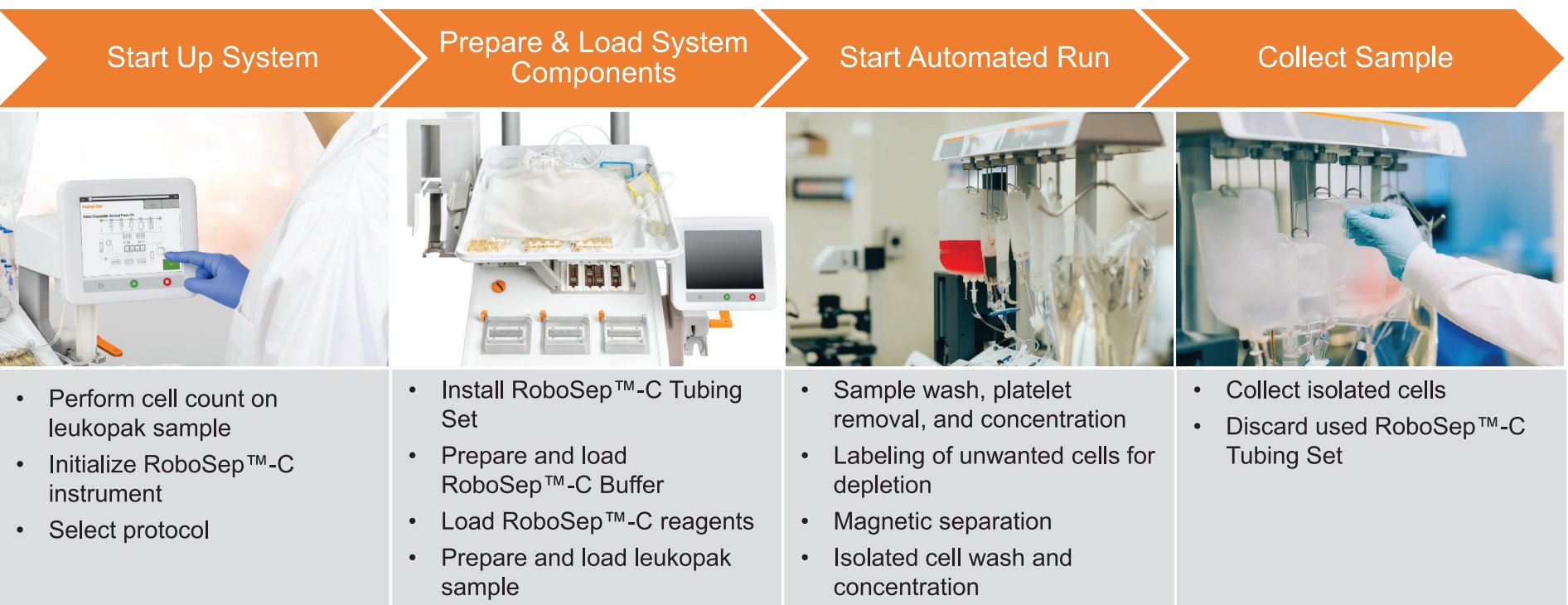


FIGURE 2. RoboSep™-C Protocol for Automated Cell Isolation.

Overview of instrument setup and automated protocol for isolation of target immune cells by negative selection. To initiate a run, all system components including the RoboSep™-C Tubing Set, Buffer, Cell Isolation reagents, and the starting leukopak (2.5 - 20 x 10⁹ nucleated cells) are loaded onto the instrument. RoboSep™-C automates the cell washing and isolation protocol within the closed, sterile, single-use tubing set. Cell concentration, buffer exchange, and platelet removal are performed within the Cell Wash Cartridge of the tubing set. RoboSep™-C negative selection reagents label unwanted cells including RBCs and platelets for depletion. Estimated total protocol time is 75 minutes (25 minutes of hands-on time and 50 minutes for the automated run).

RESULTS

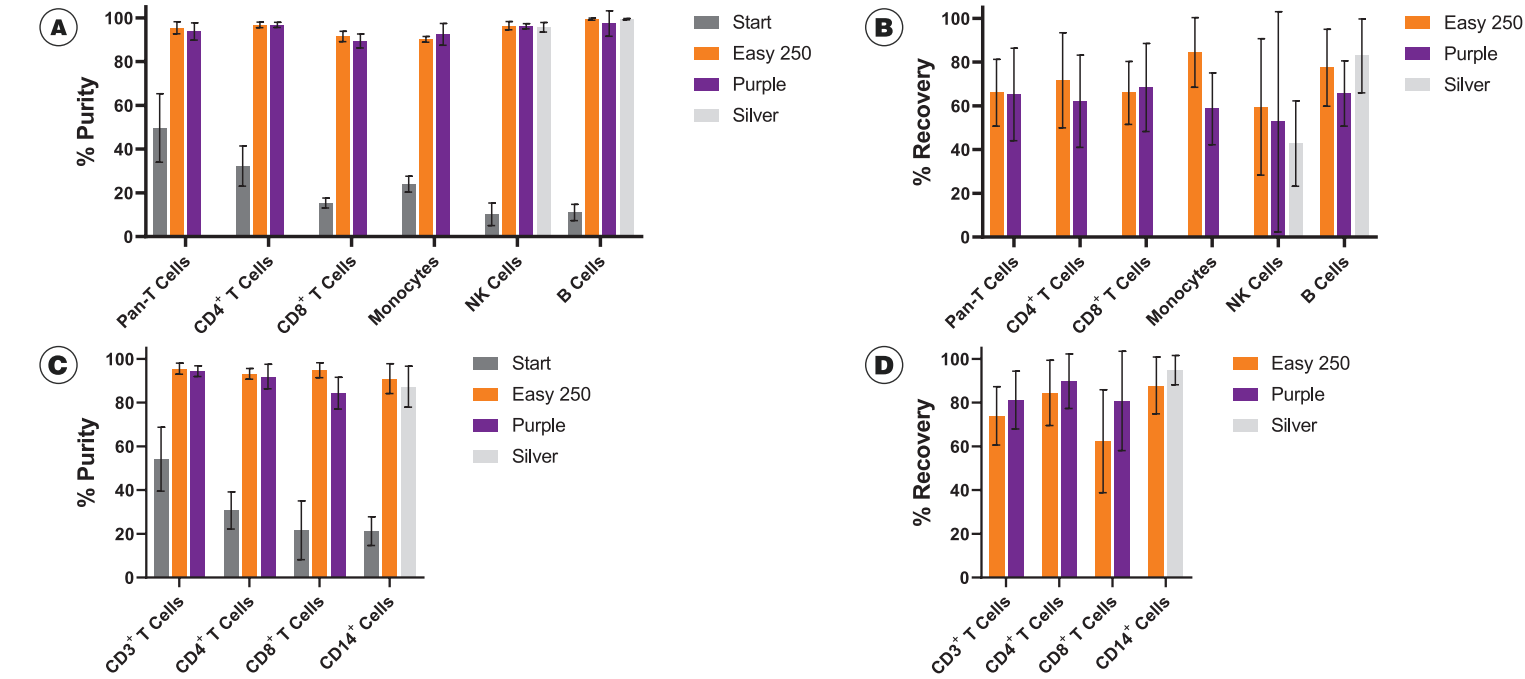


FIGURE 3. Isolation of High Purity Immune Cells from Large-Volume Samples Using the Easy 250 EasySep™ Magnet.

Immune cell subsets were isolated from processed leukopaks (washed only or washed and lysed) using various **(A - B)** negative selection or **(C - D)** positive selection EasySep™ kits. Isolations were performed using the Easy 250 EasySep™ magnet and compared to established small volume protocols using the EasySep™ magnet (purple) or “The Big Easy” EasySep™ Magnet (silver). **(A, C)** Purities of T cells (CD3⁺), CD4⁺ T cells (CD4⁺CD3⁺), CD8⁺ T cells (CD8⁺CD3⁺), monocytes (CD14⁺), NK cells (CD56⁺CD3⁺), or B cells (CD20⁺CD19⁺) within the viable CD45⁺ population were assessed by flow cytometry. **(B, D)** Recoveries are shown as the number of target cells in the isolated fraction divided by the number of target cells in the start sample x 100%. Data are shown as mean ± SD; n = 4 – 12 for each cell isolation kit.

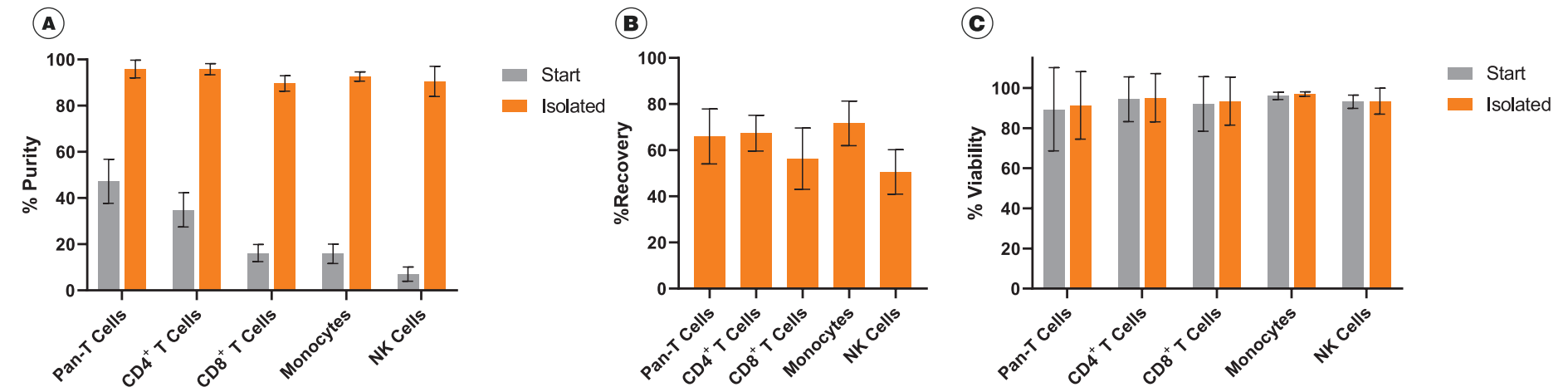


FIGURE 4. RoboSep™-C Enables Large-Scale Automated Negative Selection of Immune Cells with High Purity, Recovery, and Viability.

Pan-T cells, CD4⁺ T cells, CD8⁺ T cells, monocytes, or NK cells were isolated from fresh human peripheral blood leukopaks (2.5 - 20 x 10⁹ start cells) using the RoboSep™-C system. **(A)** Purities of T cells (CD3⁺), CD4⁺ T cells (CD4⁺CD3⁺), CD8⁺ T cells (CD8⁺CD3⁺), monocytes (CD14⁺CD16⁺), or NK cells (CD56⁺CD3⁺) within the viable CD45⁺ population were assessed by flow cytometry. **(B)** Recoveries are shown as the number of target cells in the isolated fraction divided by the number of target cells in the start sample x 100%. **(C)** Start leukopak and post-isolation viability were assessed by staining with the cell viability dye DRAQ7™ and analyzed by flow cytometry. Data are shown as mean ± SD; n = 8 – 20 for each cell isolation kit.

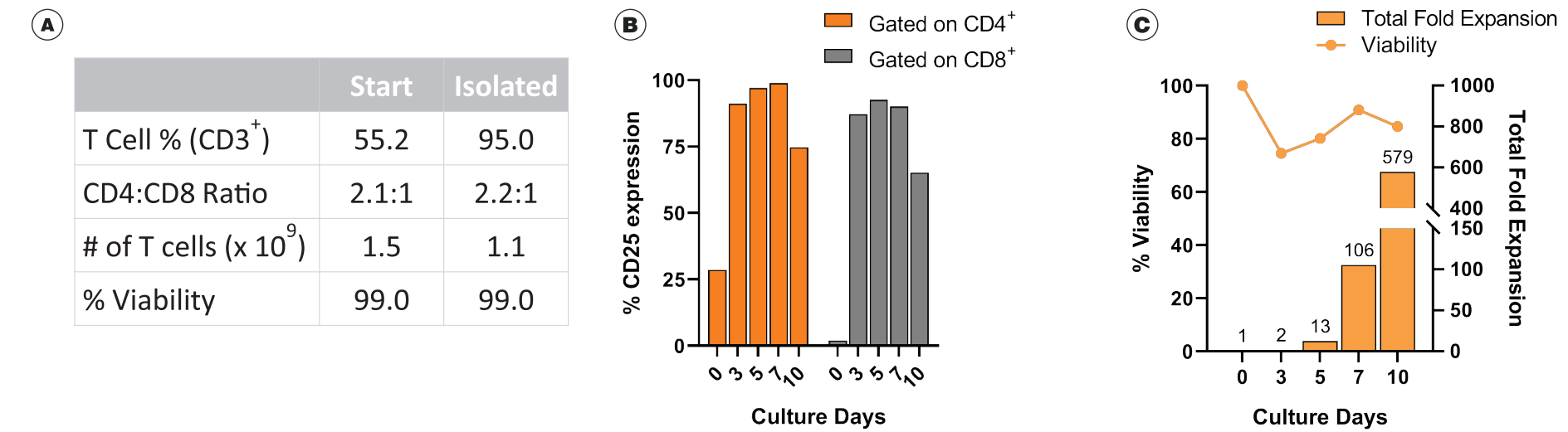


FIGURE 5. RoboSep™-C-Isolated T Cells Show Robust Activation and Expansion when Stimulated with ImmunoCult™ Human T Cell Activator in ImmunoCult™-XF T Cell Expansion Medium.

T cells were isolated from fresh human peripheral blood leukopak (2.7 x 10⁹ start cells). The isolated cells were stimulated with ImmunoCult™ Human CD3/CD28/CD2 T Cell Activator and cultured for 10 days in ImmunoCult™-XF T Cell Expansion Medium. Cells were subcultured at days 3, 5, and 7. **(A)** Analysis of start and isolated cell samples. **(B)** Expression of the activation marker CD25 was assessed by flow cytometry and **(C)** total T cell fold expansion and viability were assessed using a NucleoCounter® NC-250™ over the 10-day expansion assay. Results from a single representative experiment are shown.

Summary

- The Easy 250 EasySep™ Magnet and RoboSep™-C enable large-volume isolation of immune cells with high purity and recovery
- Isolations using the Easy 250 EasySep™ Magnet can be completed in less than 30 minutes, while the RoboSep™-C automates this procedure along with the upstream and downstream cell washing steps in a 50-minute protocol
- T cells isolated with RoboSep™-C are functional and show robust activation and expansion upon stimulation