



**THIS PRODUCT INFORMATION SHEET IS PROVIDED FOR USE WITH ROBOSEP® (SECTION A), THE PURPLE EASYSEP® MAGNET (SECTION B) OR "THE BIG EASY" SILVER EASYSEP® MAGNET (SECTION C).**

**A) FULLY AUTOMATED PROTOCOL USING ROBOSEP® (CATALOG #20000).**

This procedure is used for processing **500 µL - 8.5 mL** of sample (up to  $4.25 \times 10^8$  cells).

1. Prepare mononuclear cell suspension at a concentration of  $5 \times 10^7$  cells/mL in RoboSep® Buffer (Catalog #20104). Cells must be placed in a 14 mL (17 x 100 mm) polystyrene tube to properly fit into the RoboSep® carousel.

*Falcon™ 14 mL Polystyrene Round-Bottom Tubes (BD Biosciences, Catalog #352057) are recommended.*

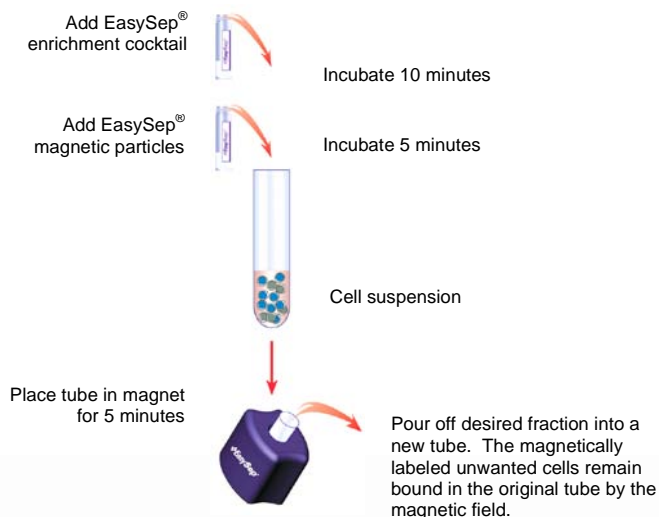
2. Select the appropriate RoboSep® protocol:
  - "Human B Cell Negative Selection 19054HLA".

If a modified RoboSep® protocol is required, please contact *STEMCELL Technologies' Technical Support* at [techsupport@stemcell.com](mailto:techsupport@stemcell.com).

3. Load the RoboSep® carousel as directed by the on-screen prompts. **Vortex EasySep® D Magnetic Particles for 30 seconds before loading. Ensure that they are in a uniform suspension with no visible aggregates.** When all desired quadrants are loaded, press the green "Run" button. All cell labeling and separation steps will be performed by RoboSep®.

4. When cell separation is complete, remove the tube containing the enriched cells from the RoboSep® carousel. Collect the enriched cells in the 50 mL tube located to the left of the tip rack. The enriched cells are now ready for use.

**MANUAL EASYSEP® PROTOCOL DIAGRAM**



**B) MANUAL EASYSEP® PROTOCOL USING PURPLE EASYSEP® MAGNET (CATALOG #18000).**

This procedure is used for processing **500 µL - 2 mL** of sample (up to  $1 \times 10^8$  cells).

1. Prepare mononuclear cell suspension at a concentration of  $5 \times 10^7$  cells/mL in recommended medium (see Notes and Tips, reverse side). Cells must be placed in a 5 mL (12 x 75 mm) polystyrene tube to properly fit into the Purple EasySep® Magnet.

*Falcon™ 5 mL Polystyrene Round-Bottom Tubes (BD Biosciences, Catalog #352058) are recommended.*

2. Add EasySep® HLA B Cell Enrichment Cocktail at **50 µL/mL cells** (e.g. for 2 mL of cells, add 100 µL of cocktail). Mix well and incubate at room temperature (15 – 25°C) for **10 minutes**.

3. Vortex EasySep® D Magnetic Particles for 30 seconds. Ensure that they are in a uniform suspension with no visible aggregates.

4. Add the EasySep® D Magnetic Particles at **75 µL/mL cells** (e.g. for 2 mL of cells, add 150 µL of particles). Mix well and incubate at room temperature (15 – 25°C) for **5 minutes**.

5. Bring the cell suspension up to a **total volume** of 2.5 mL by adding recommended medium. Mix the cells in the tube by gently pipetting up and down 2 - 3 times. Place the tube (without cap) into the magnet. Set aside for **5 minutes**.

6. Pick up the EasySep® Magnet, and in one continuous motion invert the magnet and tube, pouring off the desired fraction into a new 5 mL polystyrene tube. The magnetically labeled unwanted cells will remain bound inside the original tube, held by the magnetic field of the EasySep® magnet. Leave the magnet and the tube in inverted position for 2 - 3 seconds, then return to upright position. *Do not shake or blot off any drops that may remain hanging from the mouth of the tube.* The negatively selected, enriched cells in the new tube, are now ready for use.

**C) MANUAL EASYSEP® PROTOCOL USING "THE BIG EASY" SILVER EASYSEP® MAGNET (CATALOG #18001).**

This procedure is used for processing **500 µL - 8.5 mL** of sample (up to  $4.25 \times 10^8$  cells).

1. Prepare mononuclear cell suspension at a concentration of  $5 \times 10^7$  cells/mL in recommended medium (see Notes and Tips, reverse side). Cells must be placed in a 14 mL (17 x 100 mm) polystyrene tube to properly fit into the Silver EasySep® magnet.

*Falcon™ 14 mL Polystyrene Round-Bottom Tubes (BD Biosciences, Catalog #352057) are recommended.*

2. Add EasySep® HLA B Cell Enrichment Cocktail at **50 µL/mL cells** (e.g. for 2 mL of cells, add 100 µL of cocktail). Mix well and incubate at room temperature (15 – 25°C) for **10 minutes**.

3. Vortex EasySep® D Magnetic Particles to ensure that they are in a uniform suspension. Add the EasySep® D Magnetic Particles at **75 µL/mL cells** (e.g. for 2 mL of cells, add 150 µL of particles). Mix well and incubate at room temperature (15 – 25°C) for **5 minutes**.

4. Bring the cell suspension to a **total volume** of 5.0 mL (for  $< 2 \times 10^8$  cells) or 10 mL (for  $2 - 4.25 \times 10^8$  cells) by adding recommended medium. Mix the cells in the tube by gently pipetting up and down 2 - 3 times. Place the tube (without cap) into the magnet. Set aside for **5 minutes** at room temperature (15 – 25°C).

5. Pick up the EasySep® Magnet, and in one continuous motion invert the magnet and tube, pouring off the desired fraction into a new 14 mL polystyrene tube. The magnetically labeled unwanted cells will remain bound inside the original tube, held by the magnetic field of the EasySep® magnet. Leave the magnet and the tube inverted for 2 - 3 seconds, then return to upright position. *Do not shake or blot off any drops that may remain hanging from the mouth of the tube.* The negatively selected enriched cells in the new tube are now ready for use.

## Components:

- EasySep<sup>®</sup> HLA B Cell Enrichment Cocktail 1.0 mL
- EasySep<sup>®</sup> D Magnetic Particles 2 x 1.0 mL



NEGATIVE SELECTION

**REQUIRED EQUIPMENT:**

EasySep<sup>®</sup> Magnet (Catalog #18000), or "The Big Easy" EasySep<sup>®</sup> Magnet (Catalog #18001), or RoboSep<sup>®</sup> (Catalog #20000).

**PRODUCT DESCRIPTION AND APPLICATIONS:**

EasySep<sup>®</sup> HLA B Cell Enrichment Cocktail and EasySep<sup>®</sup> D Magnetic Particles label non-B cells for magnetic separation. These reagents are designed to enrich B cells from normal fresh or previously frozen peripheral blood mononuclear cells by depletion of non-B cells. For isolation of B cells from blood or other tissues of patients with B cell leukemia or lymphoma, or with other disease states in which B cells may express CD43, we recommend the EasySep<sup>®</sup> Human B Cell Enrichment Kit without CD43 Depletion (Catalog #19154).

**EASYSEP<sup>®</sup> LABELING OF HUMAN CELLS:**

Unwanted cells are specifically labeled with dextran-coated magnetic particles using bispecific Tetrameric Antibody Complexes (TAC). These complexes recognize both dextran and the cell surface antigen expressed on the unwanted cells (Figure 1). The small size of the magnetic dextran iron particles allows for efficient binding to the TAC-labeled cells. Magnetically labeled cells are then separated from unlabeled target cells using the EasySep<sup>®</sup> procedure (reverse side).

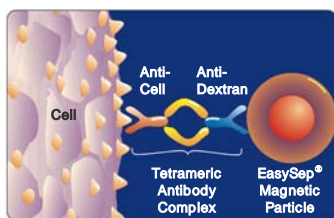


Figure 1.  
Schematic Drawing of EasySep<sup>®</sup> TAC  
Magnetic Labeling of Human Cells.

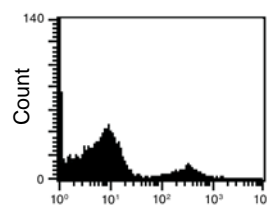
**NOTES AND TIPS:**

**PREPARING A MONONUCLEAR CELL SUSPENSION** Prepare a mononuclear cell suspension from whole peripheral blood by Ficoll-Paque<sup>™</sup> PLUS density separation (Catalog #07957). For previously frozen mononuclear cells, we recommend incubating the cells with DNase I (Catalog #07900) at a concentration of 100 µg/mL for at least 15 minutes at room temperature (15 – 25°C) prior to labeling and separation. Filter clumpy suspensions through a 30 µm mesh nylon strainer for optimal results.

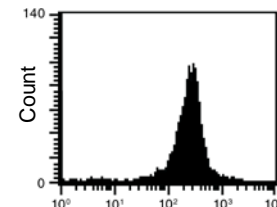
**OPTIMAL CELL NUMBER** The use of fewer than  $5 \times 10^7$  cells per separation may result in sub-optimal performance.

**RECOMMENDED MEDIUM** The recommended medium is RoboSep<sup>®</sup> Buffer (Catalog #20104), or Phosphate Buffered Saline (PBS) + 2% FBS (Catalog #07905) with 1 mM EDTA added. Medium should be Ca<sup>++</sup> and Mg<sup>++</sup> free.

**ASSESSING PURITY** Purity of B cells can be measured by flow cytometry after staining with a fluorochrome-conjugated antibody against CD19 (e.g. FITC anti-CD19, Catalog #10409), CD20 or other B cell marker.

**TYPICAL EASYSEP<sup>®</sup> HLA B CELL ENRICHMENT PROFILE:**Start: 8% CD19<sup>+</sup> CellsEnriched: 99% CD19<sup>+</sup> Cells

CD19 PE



CD19 PE

Starting with mononuclear cells, the CD19<sup>+</sup> cell content of the enriched fraction typically ranges from 95 - 99%.

**COMPONENT DESCRIPTIONS:****EASYSEP<sup>®</sup> HLA B CELL ENRICHMENT COCKTAIL****CODE #19054HC.1**

This cocktail contains a combination of monoclonal antibodies purified from hybridoma culture supernatant by affinity chromatography using Protein A or Protein G Sepharose. These antibodies are bound in bispecific Tetrameric Antibody Complexes (TAC) which are directed against cell surface antigens on human blood cells (CD2, CD3, CD14, CD16, CD36, CD43, CD56, CD66b, glycoprotein A) and dextran. The mouse monoclonal antibody subclass is IgG<sub>1</sub>. It should be noted that this product is a biological reagent, and as such cannot be completely characterized or quantified. Some variability is unavoidable.

**EASYSEP<sup>®</sup> D MAGNETIC PARTICLES****CODE #19250**

A suspension of magnetic dextran iron particles in TRIS buffer.

**STABILITY AND STORAGE:****EASYSEP<sup>®</sup> HLA B CELL ENRICHMENT COCKTAIL**

Product stable at 2 - 8°C until expiry date as indicated on label. Contents have been sterility tested. Do not freeze this product. This product may be shipped at room temperature (15 - 25°C), and should be refrigerated upon receipt.

**EASYSEP<sup>®</sup> D MAGNETIC PARTICLES**

Product stable at 2 - 8°C until expiry date as indicated on label. Contents have been sterility tested. Do not freeze this product. This product may be shipped at room temperature (15 - 25°C), and should be refrigerated upon receipt.

Ficoll<sup>™</sup> and Ficoll-Paque<sup>™</sup> PLUS are trademarks of GE Healthcare Ltd.