

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 a 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.

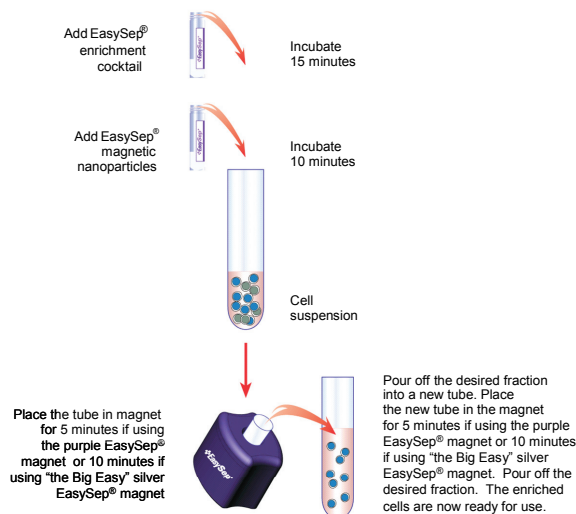
*Falcon™ 14 mL Polystyrene Round-Bottom Tubes (BD, Catalog #352057) are recommended.*

2. Select the appropriate RoboSep® protocol:
  - "Human CD4<sup>+</sup>CD127<sup>low</sup> T Cell Negative Selection 19231-high purity".

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. Mix EasySep® Magnetic Nanoparticles before loading to ensure that they are in a uniform suspension by vigorously pipetting up and down more than 5 times. Vortexing is not recommended. 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, collect the enriched cells in the 14 mL tube located to the left of the magnet in the second quadrant. 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 **250 µL - 2 mL** of sample (up to  $1 \times 10^8$  cells).

1. Prepare a mononuclear cell suspension at a concentration of  $5 \times 10^7$  cells/mL in the 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, Catalog #352058) are recommended.*

2. Add EasySep® Human CD4<sup>+</sup>CD127<sup>low</sup> T 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 **15 minutes**.
3. Mix EasySep® Magnetic Nanoparticles to ensure that they are in a uniform suspension by vigorously pipetting up and down more than 5 times. Vortexing is not recommended. Add the nanoparticles at **50 µL/mL cells** (e.g. for 2 mL of cells, add 100 µL of nanoparticles). Mix well and incubate at room temperature (15 – 25°C) for **10 minutes**.
4. Bring the cell suspension up to a **total volume** of 2.5 mL by adding the 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**.
5. 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.
6. Remove the first tube from the EasySep® Magnet and place the new tube containing the desired cells into the magnet. Set aside for **5 minutes**. 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 a mononuclear cell suspension at a concentration of  $5 \times 10^7$  cells/mL in the 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, Catalog #352057) are recommended.*

2. Add EasySep® Human CD4<sup>+</sup>CD127<sup>low</sup> T 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 **15 minutes**.
3. Mix EasySep® Magnetic Nanoparticles to ensure that they are in a uniform suspension by pipetting up and down vigorously more than 5 times. Vortexing is not recommended. Add the particles at **50 µL/ mL cells** (e.g. for 2 mL of cells, add 100 µL of nanoparticles). Mix well and incubate at room temperature (15 – 25°C) for **10 minutes**.
4. Bring the cell suspension to a **total volume** of **5.0 mL** (for  $<10^8$  cells) or **10 mL** (for  $1 - 4.25 \times 10^8$  cells) by adding the 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 **10 minutes**.
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 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 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.*
6. Remove the first tube from the EasySep® Magnet and place the new tube containing the desired cells into the magnet. Set aside for **10 minutes**. The negatively selected, enriched cells in the new tube are now ready for use.

## Components:

- EasySep<sup>®</sup> Human CD4<sup>+</sup>CD127<sup>low</sup> T Cell Enrichment Cocktail 1.0 mL
- EasySep<sup>®</sup> Magnetic Nanoparticles 2 x 1.0 mL



NEGATIVE SELECTION

**REQUIRED EQUIPMENT:**

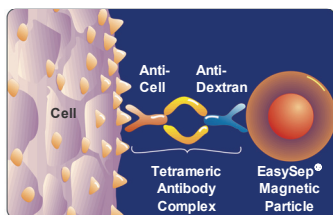
EasySep<sup>®</sup> Magnet (Catalog #18000), or "The Big Easy" EasySep<sup>®</sup> Magnet (Catalog #18001), or RoboSep<sup>®</sup> (Catalog #20000).

**PRODUCT DESCRIPTION AND APPLICATIONS:**

EasySep<sup>®</sup> Negative Selection Human CD4<sup>+</sup>CD127<sup>low</sup> T Cell Enrichment Cocktail and EasySep<sup>®</sup> Magnetic Nanoparticles label non-CD4<sup>+</sup> T cells and CD127<sup>high</sup> cells for magnetic separation. These reagents are designed to enrich CD4<sup>+</sup>CD127<sup>low</sup> T cells from fresh or previously frozen peripheral blood mononuclear cells by depletion of non-CD4<sup>+</sup> T cells and CD127<sup>high</sup> cells.

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

Unwanted cells are specifically labeled with dextran-coated magnetic nanoparticles 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.

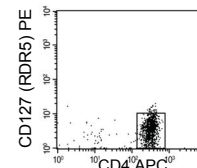
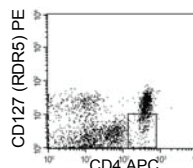
**NOTES AND TIPS:**

**PREPARING A MONONUCLEAR CELL SUSPENSION.** Prepare a mononuclear cell suspension from whole peripheral blood by Ficoll-Paque™ 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 2.5 x 10<sup>7</sup> 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). Medium should be Ca<sup>++</sup> and Mg<sup>++</sup> free.

**ASSESSING PURITY.** Purity of CD4<sup>+</sup>CD127<sup>low</sup> T cells can be measured by flow cytometry after staining with fluorochrome-conjugated anti-CD4 (e.g. FITC anti-CD4, Catalog #10403), and anti-CD127 antibodies. For CD127 staining, we recommend clone eBioRDR5 since it is not blocked by the anti-CD127 TAC present in the selection cocktail (eBioscience, Catalog #1278).

**TYPICAL EASYSEP<sup>®</sup> CD4<sup>+</sup>CD127<sup>low</sup> T CELL ENRICHMENT PROFILE:**Start: 6.9% CD3<sup>+</sup>CD4<sup>+</sup>CD127<sup>low</sup> cellsEnriched: 92% CD3<sup>+</sup>CD4<sup>+</sup>CD127<sup>low</sup> cells

Starting with peripheral blood mononuclear cells, the CD4<sup>+</sup>CD127<sup>low</sup> cell content of the enriched fraction typically ranges from 86 - 97%.

**COMPONENT DESCRIPTIONS:****EASYSEP<sup>®</sup> HUMAN CD4<sup>+</sup>CD127<sup>low</sup> T CELL ENRICHMENT COCKTAIL****CODE #19231C**

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 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> MAGNETIC NANOPARTICLES****CODE #19150.1**

A suspension of magnetic dextran iron particles in water.

**STABILITY AND STORAGE:****EASYSEP<sup>®</sup> HUMAN CD4<sup>+</sup>CD127<sup>low</sup> T 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> MAGNETIC NANOPARTICLES**

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™ and Ficoll-Paque™ PLUS are trademarks of GE Healthcare Ltd.

THIS PRODUCT IS FOR RESEARCH USE ONLY.

NOT FOR HUMAN OR ANIMAL DIAGNOSTIC, THERAPEUTIC OR COMMERCIAL USE.

WWW.STEMCELL.COM

IN NORTH AMERICA TOLL-FREE T. 1 800 667 0322 TOLL-FREE F. 1 800 567 2899 T. 1 604 877 0713 F. 1 604 877 0704 E. INFO@STEMCELL.COM

IN EUROPE T. +33 (0)4 76 04 75 30 F. +33 (0)4 76 18 99 63 E. INFO.EU@STEMCELL.COM

IN AUSTRALIA TOLL-FREE T./F. 1 800 060 350 T. 07 5474 5042 E. INFO.AUS@STEMCELL.COM

V1.1.0  
AUGUST 2009 | #29148
 Printed on recycled paper.