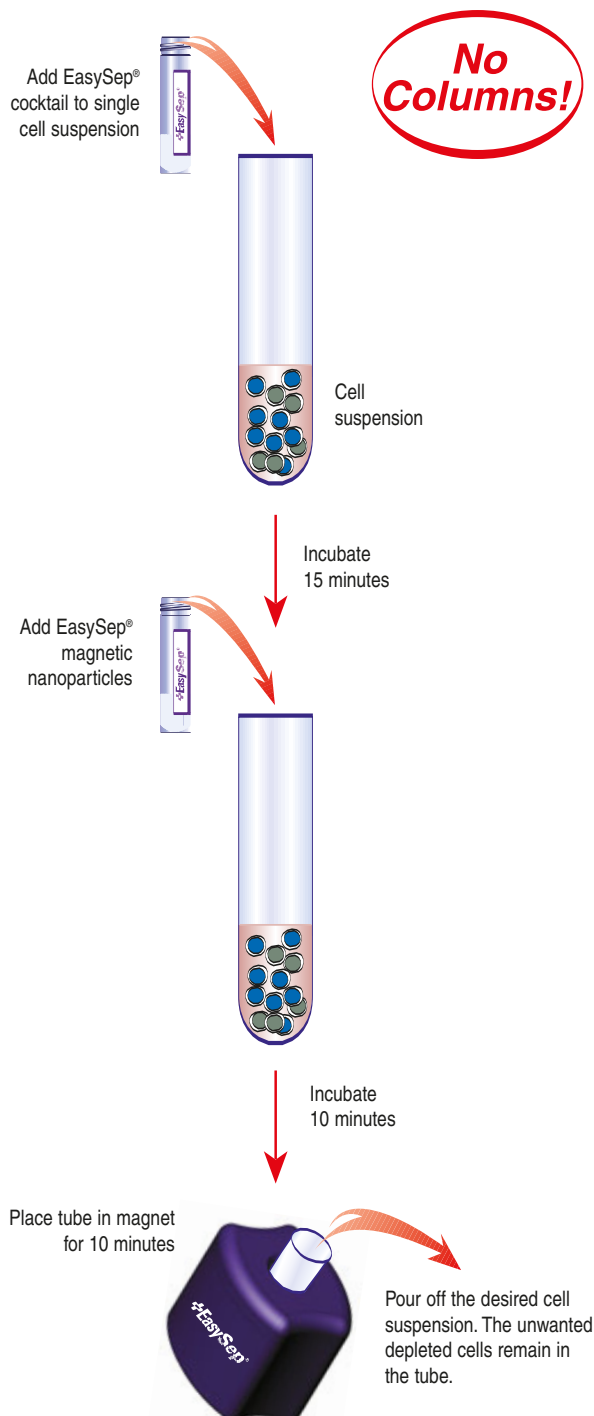


## +EasySep<sup>®</sup> Procedure:

**Note:** This procedure is designed for use with the EasySep<sup>®</sup> Magnet (18000). If using "The Big Easy" EasySep<sup>®</sup> Magnet (18001), please refer to [www.stemcell.com/technical/18001-PIS.pdf](http://www.stemcell.com/technical/18001-PIS.pdf) for additional instructions.



PROCEDURE

**Positive Selection**

# +EasySep<sup>®</sup>

## Human Cell Depletion Procedure

1. Prepare nucleated cell suspension at a concentration of  $1.25 \times 10^8$  cells/mL in recommended medium (See Notes and Tips). Cells must be placed in a 12 x 75 mm polystyrene tube to properly fit into the EasySep<sup>®</sup> Magnet. **Do not exceed a volume of 2.0 mL (i.e.  $2.5 \times 10^8$  cells) per tube.** For samples containing  $1.25 \times 10^7$  cells or fewer, resuspend in 100  $\mu$ L.  
*Falcon<sup>®</sup> 5 mL Polystyrene Round-Bottom Tubes (Becton Dickinson, Catalog #352058) are recommended.*
2. Add EasySep<sup>®</sup> Positive Selection Cocktail at 250  $\mu$ L/mL cells (eg. for 1 mL of cells add 250  $\mu$ L of cocktail). Mix well and incubate at room temperature for 15 minutes.
3. Mix EasySep<sup>®</sup> Magnetic Nanoparticles by pipetting up and down vigorously more than 5 times to ensure that they are in a uniform suspension. Vortexing is not recommended. Add the particles at 125  $\mu$ L/mL cells (eg. for 1 mL of cells add 125  $\mu$ L of nanoparticles). Mix well and incubate at room temperature for 10 minutes.
4. Bring the cell suspension 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 **ten** minutes.
5. Pick up the EasySep<sup>®</sup> Magnet, and in one continuous motion invert the magnet and tube, pouring off the supernatant fraction into a new 12 x 75 mm polystyrene tube. 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. Note:* To avoid spilling, the supernatant can be pipetted off to a new 5 mL tube or poured off into a 14 mL tube and then transferred to a 5 mL tube for the second round of separation.
6. Remove the empty tube from the EasySep<sup>®</sup> Magnet and place the new tube containing the supernatant fraction into the magnet. Set aside for **ten** minutes.
7. Repeat Step 5. The depleted cell suspension is now ready for use.

### StemCell Technologies

#### 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](mailto:info@stemcell.com)  
[www.stemcell.com](http://www.stemcell.com)

#### In the United Kingdom

Tel: +44.(0).20.7537.7565  
 Fax: +44.(0).20.7515.5408  
 Toll Free within United Kingdom:  
 Tel: 0800.731.27.14  
 Fax: 0800.731.27.13  
 e-mail: [info@stemcellgb.com](mailto:info@stemcellgb.com)  
[www.stemcellgb.com](http://www.stemcellgb.com)

#### In Europe

Tel: +33.(0).4.76.04.75.30  
 Fax: +33.(0).4.76.18.99.63  
 e-mail: [info@stemcellfrance.com](mailto:info@stemcellfrance.com)

Version 2.1  
 July 2004

**FOR RESEARCH USE ONLY**

**#28898**