

DIRECTIONS FOR USE:

Follow the instructions provided with the EasySep[®] cocktail of your choice.

The EasySep[®] Magnet is designed to hold a 12 x 75 mm polystyrene tube (we recommend Falcon[®] 5 mL Tubes, Becton Dickinson Catalog# 352058).

The size of the o-ring was selected for polystyrene tubes. Polystyrene tubes were chosen because recovery is typically better than with polypropylene tubes. Some customers may prefer to use polypropylene tubes because monocytes may adhere to polystyrene. If polypropylene tubes are used, we advise that the diameter of the opening be reduced by inserting a small piece of rubber band between the o-ring and the housing.

CAUTION:

The EasySep[®] Magnet generates a strong magnetic field. Keep away from pacemakers, magnets, computer disks, watches, and other objects that respond to magnetic fields.

PRODUCT DESCRIPTION:

A sealed anodized aluminum body housing rare earth magnetic elements.

CARE OF MAGNET:

After completing a cell separation, wipe the magnet clean with distilled water and then with alcohol to dry the magnet. Do not clean with bleach.

PROCEDURE

Magnets

+EasySep[®]



EasySep[®]
Magnet

CATALOG #18000

FOR RESEARCH USE ONLY

In North America

Tel: 1.604.877.0713
Fax: 1.604.877.0704
NA Toll Free Tel: 1.800.667.0322
NA Toll Free Fax: 1.800.567.2899
e-mail: info@stemcell.com
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
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

StemCell Technologies

Version 1.2.1
November 2004

28922