

# Lymphocyte separation media

## Instruction manual

Product	Size	Catalog number
Lymphocyte Separation Medium 1077	500 ml	C-44010

### Product description

Lymphocyte Separation Medium 1077 is intended for the separation of the portion of vital mononuclear cells from whole blood, buffy coats, bone marrow and several other starting materials, e.g., crude cell preparations, by means of low density gradient centrifugation. Lymphocyte Separation Medium has a density of 1.077 g/ml at 20°C. It is sterile, ready-to-use and based on Ficoll 400 and sodium diatrizoate providing optimal physiological parameters and low cytotoxicity.

### Instruction of use

1. Fill Lymphocyte Separation Medium 1077 in a tube. Use 3 ml for a 15 ml tube and 20 ml for a 50 ml tube.

2. Hold the tube in a 45° angle and carefully load the cell suspension to be separated on top of the separation medium.

**Note:** Perform this step slowly but constantly. Take care that the two phases do not mix and a sharp border is present.

For best results, use an electrical pipet aid with EX-function allowing the slow and constant efflux of the cell suspension from the pipet.

3. Centrifuge the tube at 440 x g for 40 minutes at room temperature without brake.

4. Carefully aspirate most of the supernatant without disturbing the layer of mononuclear cells in the interphase.

5. Aspirate the ring of mononuclear cells from the interphase using a pipet. Keep the volume as small as possible.

6. Wash the mononuclear cells with PBS/0.1% HSA or BSA and centrifuge at 360 x g for 10 minutes.

7. Wash the mononuclear cells with PBS/0.1% HSA or BSA and centrifuge at 200 x g for 10 minutes. Repeat this step.

8. The mononuclear cell preparation is ready.

### Storage and stability

Lymphocyte Separation Medium 1077 is light-sensitive. Store the medium at room temperature in the dark immediately after arrival. Do not freeze. If stored properly, the product is stable until the expiry date stated on the label.

### Quality control

All lots of PromoCell Lymphocyte Separation Medium 1077 are subjected to comprehensive quality control tests. Each lot is routinely tested for function and physical parameters such as osmolality and pH level. Approved in-house lots are used as a reference.

In addition, all lots have been tested for the absence of microbial contaminants (fungi and bacteria).

### Intended use

The products are for *in vitro* research use only and not for diagnostic or therapeutic procedures. For safety precautions please see appropriate MSDS.

**PromoCell GmbH**  
Sickingenstr. 63/65  
69126 Heidelberg  
Germany

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

**USA/Canada**

Phone: 1 - 866 - 251 - 2860 (toll free)  
Fax: 1 - 866 - 827 - 9219 (toll free)

**Deutschland**

Telefon: 0800 - 776 66 23 (gebührenfrei)  
Fax: 0800 - 100 83 06 (gebührenfrei)

**France**

Téléphone: 0800 - 90 93 32 (ligne verte)  
Téléfax: 0800 - 90 27 36 (ligne verte)

**United Kingdom**

Phone: 0800 96 03 33 (toll free)  
Fax: 0800 169 85 54 (toll free)

**Other Countries**

Phone: +49 6221 - 649 34 0  
Fax: +49 6221 - 649 34 40

© PromoCell GmbH