

# Mononuclear Cell Medium

## Instruction Manual

| Product  | Size   | Catalog Number |
|--|--------|----------------|
| Mononuclear Cell Medium (Ready-to-use) – Includes Basal Medium and SupplementMix | 500 ml | C-28030        |

### Recommended for

- Human Mononuclear Cells from Cord Blood (hMNC-CB)
- Human Mononuclear Cells from Peripheral Blood (hMNC-PB)

### Product Description

Our Mononuclear Cell Medium is a very rich short-term maintenance medium, which provides utmost cell viability for Mononuclear Cells from umbilical cord blood, adult peripheral blood and other sources, e.g. bone marrow. It is suitable to maintain Mononuclear Cells for up to 48 hours before proceeding with your experiments.

The Medium consists of a bottle of Basal Medium and one vial of SupplementMix. Adding the SupplementMix to the Basal Medium results in the complete Medium.

### Supplementation Details

The Mononuclear Cell Medium contains all the growth factors and supplements necessary for optimal short-term maintenance of

mononuclear cells. The Mononuclear Cell Medium does not contain antibiotics or antimycotics and is formulated for use in an incubator with an atmosphere of 5% CO<sub>2</sub>.

### Preparation of the Supplemented Medium for Use

Thaw the supplements at 15–25°C. Aseptically mix the supplement solutions by carefully pipetting up and down. Then, transfer the entire content of each supplement to the Basal Medium. Close the bottle and swirl gently until a homogenous mixture is formed.

### Storage and Stability

Store the Basal Medium at 4–8°C in the dark and the SupplementMix at -20°C immediately after arrival. Do not freeze the Basal Medium. If stored properly, the products are stable until the expiry date stated on the label. After adding the supplements to the Basal Medium, the shelf life of the complete medium is 6 weeks at 4–8°C. Do not freeze the complete medium. For use, pre-warm only an aliquot of the complete medium at

15–25°C and keep the remaining medium refrigerated at 4–8°C.

**Note:** The SupplementMix is delivered thawed and can be frozen after arrival without losing any activity.

### Quality Control

All lots of PromoCell Mononuclear Cell Medium are subjected to comprehensive quality control tests using primary human mononuclear cells from cord blood and adult peripheral blood. Each lot is checked for unaltered short-term maintenance of highly viable mononuclear cells. Approved in-house lots of media are used as a reference. In addition, all lots of media have been tested for the absence of microbial contaminants (fungi, bacteria, mycoplasma).

### Intended Use

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

---

If you require special media modifications, we offer a custom media service starting at 10 bottles per order.  
Contact us at [info@promocell.com](mailto:info@promocell.com) to find out more.

---

**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