

# Hematopoietic Progenitor Cell Expansion Media XF

## Instruction manual

### Hematopoietic Progenitor Cell Expansion Media XF

Product	Size	Catalog number
Hematopoietic Progenitor Cell Expansion Medium XF	500 ml	C-28021
Hematopoietic Progenitor Cell Expansion Medium XF, phenol red-free	500 ml	C-28022

### Cytokine Mix E for HPC Expansion Media XF

Product	Size	Catalog number
Cytokine Mix for HPC Expansion Media XF	1 ml (sufficient for 100 ml Medium)	C-39890
	5 ml (sufficient for 500 ml Medium)	C-39891

#### Recommended for

- Human Mononuclear Cells (hMNC)
- Human CD34<sup>+</sup> Progenitor Cells (hCD34<sup>+</sup>-CB)
- Human Hematopoietic Progenitor Cells (HPCs)

#### Product description

Our Hematopoietic Progenitor Cell (HPC) Expansion Media XF provide a highly optimized, serum-free and xeno-free culture system allowing for superior expansion performance of human hematopoietic progenitor cells, e.g., CD34<sup>+</sup> and CD133<sup>+</sup> cells. The HPC Expansion Media XF combined with Cytokine Mix E typically promotes over 300-fold expansion of the total cell population. Depending on the donor, expansions exceeding 2000-fold can be achieved. After 2 weeks of expansion about 20–30% of the population express CD34<sup>+</sup>, indicating a 50–200 fold expansion of CD34<sup>+</sup> progenitor cells.

Cytokine Mix E is a ready-to-use mixture of growth factors optionally available for the

HPC Expansion Media XF. The serum-free and xeno-free formula contains recombinant human TPO, SCF, FLT-3 ligand and IL-3.

With traditional serum-containing media, the majority of HPC lose their functionality during the expansion process. Our optimized serum-free and xeno-free formulation allows for an expanded cell population that is enriched with functional progenitor cells capable of CFC-formation in appropriate *in vitro* assays.

#### Supplementation details

The HPC Expansion Media XF must be supplemented with additional growth factors by the user, e.g., the optionally available Cytokine Mix E (C-39890, 1 ml or C-39891, 5 ml) containing recombinant human TPO, SCF, FLT-3 ligand and IL-3. However, the user may customize the media by supplementation with growth factors of his choice.

The HPC Expansion Media XF do not contain antibiotics or antimycotics and are 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.

For the HPC Expansion Media XF additionally thaw the Cytokine Mix E at 15–25°C. Aseptically mix the solutions by carefully pipetting up and down. Then, transfer the entire content to the referring amount of complete Medium. Close the bottle and swirl gently until a homogenous mixture is formed. One milliliter of Cytokine Mix E is sufficient for the supplementation of 100 ml HPC Expansion Medium XF. Alternatively, supplement the HPC Expansion Media XF with growth factors of your choice.

## Storage and stability

Store the basal medium at 4 to 8°C in the dark, store the SupplementMix and Cytokine Mix 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 SupplementMix to the basal medium, the shelf life of the complete medium is 6 weeks at 4 to 8°C. Medium supplemented with Cytokine Mix E should be

used within 2 weeks after addition of the cytokines. Do not freeze the complete medium. For use, pre-warm only an aliquot of the complete medium and keep the remaining medium refrigerated at 4 to 8°C.

## Quality control

Each lot of PromoCell HPC Expansion Media XF is tested for the ability to support the expansion of hematopoietic progenitor cells.

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

## Intended use

The products are 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

**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

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

© PromoCell GmbH