

## Instruction manual

Product	Size	Catalog number
Cryo-SFM Plus	30 ml 125 ml	C-29920 C-29922
Cryo-SFM Plus, phenol red-free	30 ml 125 ml	C-29930 C-29932

### Product description

Our Cryo-SFM Plus are defined, animal component-free, and protein-free cryopreservation media. The optimized formulation is based on methylcellulose, DMSO and other cryoprotectants. Cryo-SFM Plus and Cryo-SFM Plus, phenol red-free media provide excellent results with all types of cells including primary human cells, stem cells and cell lines. Cells frozen in Cryo-SFM Plus or Cryo-SFM Plus, phenol red-free exhibit superior viability, attachment, and subsequent growth performance after thawing.

### Instruction of use

Detach adherent cells using the DetachKit (C-41200, C-41210, C-41220). Resuspend the cells in room temperature Cryo-SFM Plus

(C-29920/C-29922) or Cryo-SFM Plus, phenol red-free (C-29930/C-29932) at a concentration of 0.1-10 million cells/ml. Transfer the cell suspension into cryovials, freeze them gradually (-1°C/minute) to -70°C or lower overnight and subsequently store the vials in liquid nitrogen.

**Note:** Each freeze/thaw cycle leads to a loss in proliferation potential and cell number, particularly when working with fragile primary cells. When freezing cells, use the lowest passage number possible.

### Storage and stability

Store Cryo-SFM Plus and Cryo-SFM Plus, phenol red-free at 4 to 8°C in the dark immediately after arrival. If stored properly, the product is stable until the expiry date stated on the label.

### Quality control

All lots of PromoCell Cryo-SFM Plus and Cryo-SFM Plus, phenol red-free are subjected to comprehensive quality control tests. Each lot is routinely tested for its performance with our primary Human Mesenchymal Stem Cells from Adipose Tissue (hMSC-AT).

In addition, all lots of Cryo-SFM Plus and Cryo-SFM Plus, phenol red-free 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  
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