

Instruction manual

Product	Size	Catalog number
Cryo-SFM Plus	30 ml 125 ml	C-29920 C-29922

Product description

Our Cryo-SFM Plus is a defined, animal component-free, and protein-free cryopreservation medium. The optimized formulation is based on methylcellulose, DMSO and other cryoprotectants. Cryo-SFM Plus provides excellent results with all types of cells including primary human cells, stem cells and cell lines. Cells frozen in Cryo-SFM Plus 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 cold or room temperature Cryo-SFM Plus 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. If freezing the cells, please do so at as low a passage number as possible.

Storage and stability

Store Cryo-SFM Plus 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 are subjected to comprehensive quality control tests. Each lot is routinely tested for its performance with primary human cells. Approved in-house lots are used as a reference.

In addition, all lots of Cryo-SFM Plus 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