

Instruction Manual

Product	Size	Catalog Number
Cryo-SFM	30 ml 125 ml	C-29910 C-29912

Product Description

Our Cryo-SFM is a defined, animal component-free, and protein-free cryopreservation medium. The optimized formulation is based on methylcellulose, DMSO and other cryoprotectants. Cryo-SFM provides excellent results with all types of cells including primary human cells, stem cells and cell lines. Cells frozen in Cryo-SFM 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 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 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 are subjected to comprehensive quality control tests. Each lot is routinely tested for its performance with primary human cells, for the absence of cytotoxicity, and for physical parameters such as osmolality and pH level. Approved in-house lots are used as a reference.

In addition, all lots of Cryo-SFM have been tested for the absence of microbial contaminants (fungi, 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