

## Instruction Manual

| Product                                   | Size                       | Catalog Number |
|---|----------------------------|----------------|
| NHEK.f single donor Pellet                | 1 million cells per pellet | C-14001        |
| NHEK adult, single donor Pellet           | 1 million cells per pellet | C-14002        |
| NHEK.f pooled Pellet                      | 1 million cells per pellet | C-14003        |
| NHEK adult pooled Pellet                  | 1 million cells per pellet | C-14004        |
| HFDPIC Pellet                             | 1 million cells per pellet | C-14005        |
| HUVEC Growth Medium 2 single donor Pellet | 1 million cells per pellet | C-14008        |
| HUVEC Growth Medium 2 pooled Pellet       | 1 million cells per pellet | C-14009        |
| HUVEC single donor Pellet                 | 1 million cells per pellet | C-14010        |
| HUVEC pooled Pellet                       | 1 million cells per pellet | C-14011        |
| HUVEC pre-screened Pellet                 | 1 million cells per pellet | C-14012        |
| HUAEC Pellet                              | 1 million cells per pellet | C-14013        |
| HDMEC Pellet                              | 1 million cells per pellet | C-14015        |
| HDMEC adult Pellet                        | 1 million cells per pellet | C-14016        |
| HDMEC pre-screened Pellet                 | 1 million cells per pellet | C-14017        |
| HDBEC Pellet                              | 1 million cells per pellet | C-14018        |
| HDBEC adult Pellet                        | 1 million cells per pellet | C-14019        |
| HDLEC Pellet                              | 1 million cells per pellet | C-14020        |
| HDLEC adult Pellet                        | 1 million cells per pellet | C-14021        |
| HCAEC Pellet                              | 1 million cells per pellet | C-14022        |
| HAoEC Pellet                              | 1 million cells per pellet | C-14023        |
| HPAEC Pellet                              | 1 million cells per pellet | C-14024        |
| HSaVEC Pellet                             | 1 million cells per pellet | C-14025        |
| HPMEC Pellet                              | 1 million cells per pellet | C-14027        |
| HUtMEC Pellet                             | 1 million cells per pellet | C-14028        |
| HCMEC Pellet                              | 1 million cells per pellet | C-14029        |

| Product  | Size                       | Catalog Number |
|--|----------------------------|----------------|
| NHDF Pellet                                    | 1 million cells per pellet | C-14030        |
| NHDF adult Pellet                              | 1 million cells per pellet | C-14031        |
| HPF Pellet                                     | 1 million cells per pellet | C-14035        |
| HCF Pellet                                     | 1 million cells per pellet | C-14036        |
| HAoAF Pellet                                   | 1 million cells per pellet | C-14037        |
| HUF Pellet                                     | 1 million cells per pellet | C-14038        |
| NHEM.f Pellet                                  | 1 million cells per pellet | C-14040        |
| NHEM.f M3 Pellet                               | 1 million cells per pellet | C-14044        |
| NHEM M3 adult Pellet                           | 1 million cells per pellet | C-14045        |
| NHEK.f Growth Medium 3 single donor Pellet     | 1 million cells per pellet | C-14046        |
| NHEK adult Growth Medium 3 single donor Pellet | 1 million cells per pellet | C-14047        |
| NHEK.f Growth Medium 3 pooled Pellet           | 1 million cells per pellet | C-14048        |
| NHEK adult Growth Medium 3 pooled Pellet       | 1 million cells per pellet | C-14049        |
| HUASMC Pellet                                  | 1 million cells per pellet | C-14050        |
| HPASMC Pellet                                  | 1 million cells per pellet | C-14051        |
| HCASMC Pellet                                  | 1 million cells per pellet | C-14052        |
| HAoSMC Pellet                                  | 1 million cells per pellet | C-14053        |
| HBSMC Pellet                                   | 1 million cells per pellet | C-14055        |
| HTSMC Pellet                                   | 1 million cells per pellet | C-14056        |
| HUtSMC Pellet                                  | 1 million cells per pellet | C-14059        |
| SkMC Pellet                                    | 1 million cells per pellet | C-14060        |
| HNEpC Pellet                                   | 1 million cells per pellet | C-14062        |
| HBepC Pellet                                   | 1 million cells per pellet | C-14063        |
| HTEpC Pellet                                   | 1 million cells per pellet | C-14064        |
| HSAEpC Pellet                                  | 1 million cells per pellet | C-14065        |
| HREpC Pellet                                   | 1 million cells per pellet | C-14067        |
| HRCEpC Pellet                                  | 1 million cells per pellet | C-14068        |
| HCH Pellet                                     | 1 million cells per pellet | C-14070        |
| HOB Pellet                                     | 1 million cells per pellet | C-14071        |
| HWP subcutaneous Pellet                        | 1 million cells per pellet | C-14072        |

| Product                         | Size                       | Catalog Number |
|---------------------------------|----------------------------|----------------|
| HWP visceral Pellet             | 1 million cells per pellet | C-14073        |
| HCM Pellet                      | 1 million cells per pellet | C-14080        |
| hMSC-BM Pellet                  | 1 million cells per pellet | C-14090        |
| hMSC-UC Pellet                  | 1 million cells per pellet | C-14091        |
| hMSC-AT Pellet                  | 1 million cells per pellet | C-14092        |
| hPC-PL Pellet                   | 1 million cells per pellet | C-14095        |
| hMNC-CB single donor Pellet     | 1 million cells per pellet | C-14096        |
| hMoCD14+-PB single donor Pellet | 1 million cells per pellet | C-14110        |
| hMNC-PB single donor Pellet     | 1 million cells per pellet | C-14098        |

### Product Description

Our Cell Pellets are stored in RNAlater® – an aqueous protection solution that rapidly permeates the cells and stabilizes cellular DNA, RNA, and proteins and protects them from degradation. This makes our Cell Pellets very useful for a wide range of research applications like gene expression analysis, cell characterization studies, and protein characterization experiments.\*

**Note:** RNAlater® denatures proteins. Therefore, proteins will be suitable for e.g. Western blotting but not for applications that require native proteins or viable cells. Because of the density of RNAlater® solution, greater centrifugal forces are required to pellet cells. We recommend to spin down the cells at 5000 x g at 4°C. Alternatively, if no pellet is visible after centrifugation, RNA can be purified directly from the RNAlater® solution. This can be done by adding 2 ml of 10x lysis buffer, and proceeding normally.

Each Cell Pellet is made from 1 million cells and is dissolved in 200 µl RNAlater®. The Cell Pellets are produced from released cell lots and have the same characteristics, i.e. expression profile or cell type specific antigens. On request, viable cells from the same lot are available for cultivation.

### Storage and Stability

The Cell Pellets can be stored indefinitely at -20°C, and for up to 1 month at 4°C.

**Note:** The Cell Pellets will not freeze at -20°C. Always centrifuge product briefly before opening vial.

### Quality Control

Rigid quality control tests are performed for each lot of the corresponding viable cells. They are tested for cell viability, cell morphology, growth performance and for cell-type specific markers.

In addition, all cells have been tested for the absence of HIV-1, HIV-2, HBV, HCV, HTLV-1, HTLV-2 and microbial contaminants (fungi, bacteria, and mycoplasma).

### Intended Use

Our Cell Pellets are for in vitro research use only and not for diagnostic or therapeutic procedures.

### Warning

Although tested negative for HIV-1, HIV-2, HBV, HCV, HTLV-1 and HTLV-2, the cells – like all products of human origin – should be handled as potentially infectious. No test procedure can completely guarantee the absence of infectious agents.

*Follow appropriate safety precautions!*

\* More information on RNAlater protocols and compatible kits at [www.thermofisher.com/de/de/home/brands/product-brand/rnalater.html](http://www.thermofisher.com/de/de/home/brands/product-brand/rnalater.html)

**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