

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
HFDPK 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 for 10 minutes. 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 and cell pellets from the same lot can be offered.

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/malater.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