

Instruction manual

Product	Size	Catalog number
Normal Human Dermal Fibroblasts (NHDF) juvenile foreskin	500,000 cryopreserved cells 500,000 proliferating cells	C-12300 C-12350
Normal Human Dermal Fibroblasts (NHDF) adult donor	500,000 cryopreserved cells 500,000 proliferating cells	C-12302 C-12352
Human Pulmonary Fibroblasts (HPF)	500,000 cryopreserved cells 500,000 proliferating cells	C-12360 C-12361
Human Aortic Adventitial Fibroblasts (HAoAF)	500,000 cryopreserved cells 500,000 proliferating cells	C-12380 C-12381
Human Cardiac Fibroblasts (HCF)	500,000 cryopreserved cells 500,000 proliferating cells	C-12375 C-12377
Human Uterine Fibroblasts (HUF)	500,000 cryopreserved cells 500,000 proliferating cells	C-12385 C-12386

Product description

Derived from the mesoderm, connective tissue is involved in maintaining the structural integrity of the body. The predominant cells found in connective tissue are fibroblasts, which continuously secrete various components of the extracellular matrix. Depending on the site of origin and physiological status, fibroblasts may exhibit different morphological phenotypes and various functional properties.

We offer a range of Human Fibroblasts produced at our cell culture facility. The cells are isolated from normal human juvenile foreskin or adult skin from different locations. Additionally, we offer fibroblasts from adult lung tissue and aortic adventitial tissue.

Shortly after isolation, all our Human Fibroblasts are cryopreserved using our proprietary, animal-component free, and protein-free cryopreservation medium (for details, please refer to the Certificate of Analysis).

Each cryovial contains more than 500,000 viable cells after thawing. Thawing and seeding results in passage 2.

Proliferating cell cultures are made from 500,000 cryopreserved cells that have been thawed and cultured for three days in our hands and shipped as growing cultures.

Quality control

We perform rigid quality control tests for each lot of Human Fibroblasts.

The cells are tested for cell morphology, adherence rate and cell viability. Flow cytometric analyses of the fibroblast specific antigen CD90 are performed for each lot. For tested marker details, please refer to the Certificate of Analysis. Growth performance is tested through multiple passages up to 10 or 15 population doublings (PD) under culture conditions without antibiotics or antimycotics (see page 5).

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 Human Fibroblasts 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!

After delivery, cryopreserved cells should be stored in liquid nitrogen or seeded directly (see page 2). Proliferating cells must be processed immediately (see page 3).

Protocol for cryopreserved cells

Straight after arrival, store the cryopreserved cells in liquid nitrogen, or seed them immediately.

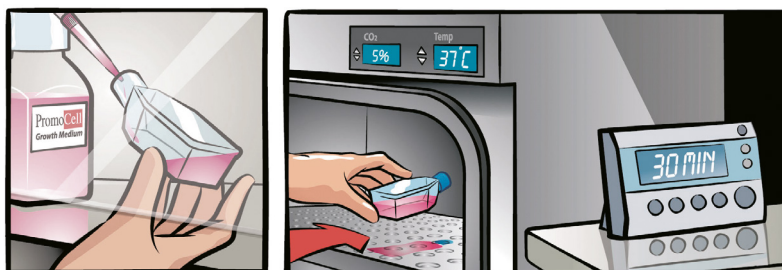
Note: Storage at -80°C is not sufficient for cell preservation and causes irreversible cell damage.

Use aseptic techniques and a laminar flow bench.

1

Prepare the medium

Calculate the required culture surface area according to the plating density (see page 5) and the lot-specific cell numbers stated on the certificate of analysis. Fill the appropriate volume of PromoCell Growth Medium (at least 9 ml per vial of cells) in cell culture vessels. Place the vessels in an incubator (37°C , 5% CO_2) for 30 minutes.



2

Thaw the cells

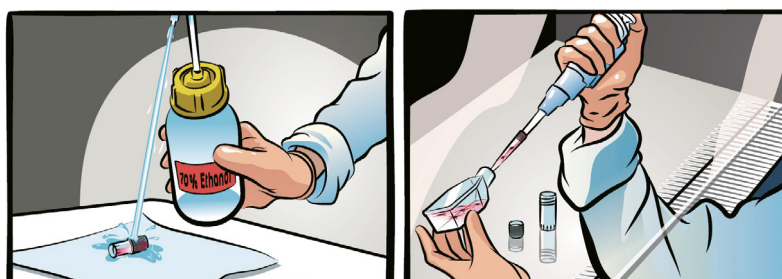
Remove the cryovial from the liquid nitrogen container and immediately place it on dry ice – even for short transportation. Under a laminar flow bench, briefly twist the cap a quarter turn to relieve pressure, then retighten. Immerse the vial in a water bath (37°C) up to the height of the screw cap for 2 minutes. Ensure that no water enters the thread of the screw cap.



3

Disinfect the vial and seed the cells

Thoroughly rinse the cryovial with 70% ethanol under a laminar flow bench. Then, aspirate the excess ethanol from the thread area of the screw cap. Open the vial and transfer the cells to a cell culture vessel containing the pre-warmed medium from step 1.



4

Incubate the cells

Place the vessel in an incubator (37°C , 5% CO_2) for cell attachment. Replace the medium after 16–24 hours and every two to three days thereafter. The cells should be subcultured, according to the subcultivation protocol (see page 4), once they have reached 70–90% confluency.



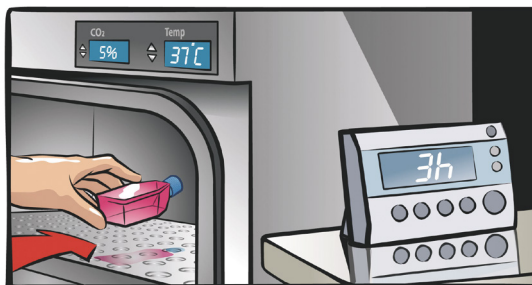
Protocol for proliferating cells

Start immediately after delivery.
Use aseptic techniques and a laminar flow bench.

1

Incubate the cells

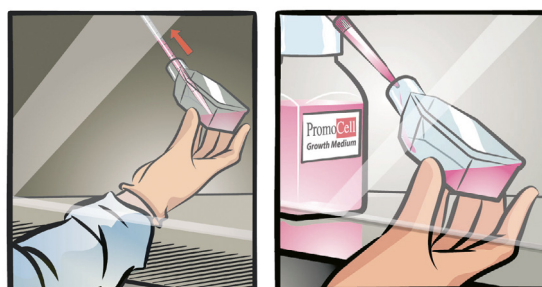
Unpack the culture vessel, do not open the cap, and immediately place it in an incubator (37°C, 5% CO₂) for 3 hours to allow the cells to recover from transportation.



2

Replace the transport medium

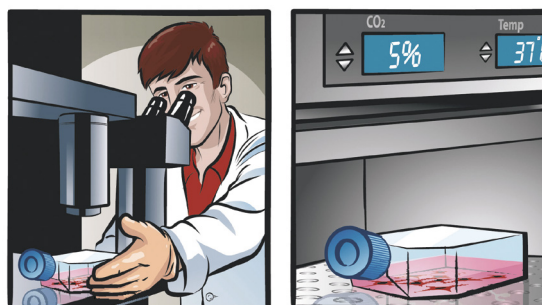
Carefully open the vessel, rinse the inner side of the cap with 70% ethanol, and let air dry. Aspirate the transport medium from the vessel. Add 10 ml of the appropriate PromoCell Cell Growth Medium.



3

Check and incubate the cells

Check the cell density. Open the cap half a turn and place the vessel in an incubator (37°C, 5% CO₂). Change the medium every two to three days. The cells should be subcultured, according to the subcultivation protocol (see page 4), once they have reached 70–90% confluency.



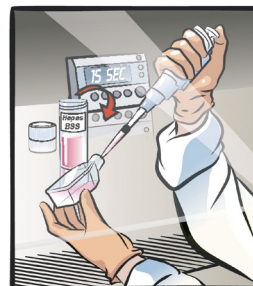
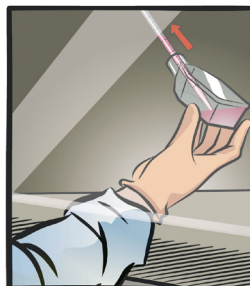
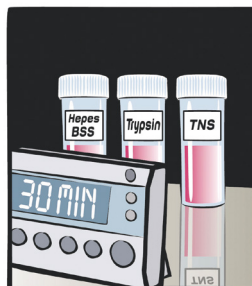
Subcultivation protocol

Use aseptic techniques and a laminar flow bench.

1

Prepare the reagents and wash the cells

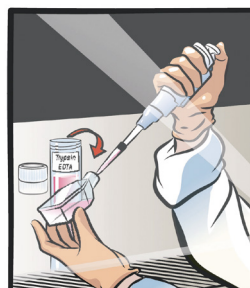
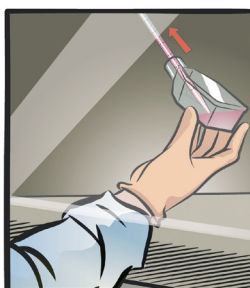
Place the PromoCell DetachKit at room temperature for at least 30 minutes to adjust the temperature of the reagents. Carefully aspirate the medium from the culture vessel. Add 100 μ l Hepes BSS Solution per cm^2 of vessel surface to wash the cells and agitate the vessel carefully for 15 seconds.



2

Detach the cells

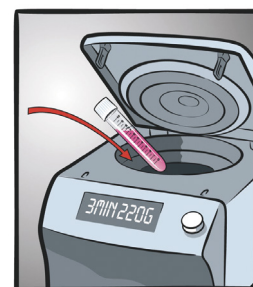
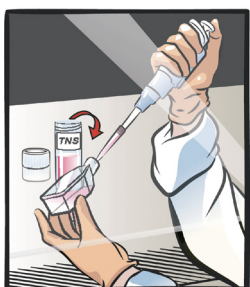
Carefully aspirate the Hepes BSS from the culture vessel. Add 100 μ l Trypsin/EDTA Solution per cm^2 of vessel surface. Note: We recommend detaching the cells at room temperature. Close the vessel and examine the cells under a microscope. When the cells start to detach, gently tap the side of the vessel to loosen the remaining cells.



3

Neutralize the trypsin and harvest the cells

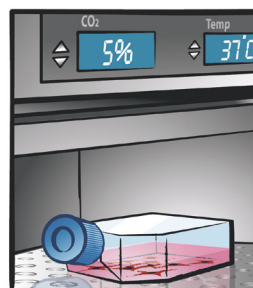
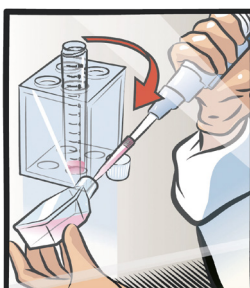
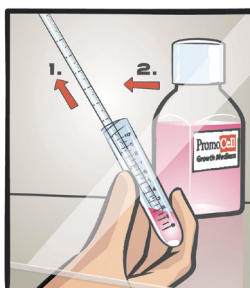
Add 100 μ l Trypsin Neutralization Solution per cm^2 of vessel surface and gently agitate. Carefully aspirate the cell suspension and transfer it to a centrifugation tube. Spin down the cells for 3 minutes at 220 x g.



4

Incubate the cells

Discard the supernatant (step 1), add 1 ml of the appropriate PromoCell Cell Growth Medium (step 2), and re-suspend the cells by carefully pipetting up and down. Plate the cells according to the recommended seeding density in new cell culture vessels containing prewarmed PromoCell Cell Growth Medium. Place the vessels in an incubator (37°C, 5% CO_2) and change the media every two to three days.



Specifications

Product	Recommended culture media*	Plating density	Passage after thawing	Marker	Population doublings
Normal Human Dermal Fibroblasts (NHDF), juvenile foreskin	C-23010	3,500 – 7,000 cells per cm ²	P2	CD90 ⁺	> 15
Normal Human Dermal Fibroblasts (NHDF), adult donor	C-23020	3,500 – 7,000 cells per cm ²	P2	CD90 ⁺	> 15
Human Pulmonary Fibroblasts (HPF)	C-23020	3,500 – 7,000 cells per cm ²	P2	CD90 ⁺	> 15
Human Aortic Adventitial Fibroblasts (HAoAF)	C-23020	3,500 – 7,000 cells per cm ²	P2	CD90 ⁺	> 10
Human Cardiac Fibroblasts (HCF)	C-23025	3,500 – 7,000 cells per cm ²	P2	CD90 ⁺ Fibronectin ⁺	> 15
Human Uterine Fibroblasts (HUF)	C-23020	3,500 – 7,000 cells per cm ²	P2	CD90 ⁺	> 15

**The catalog numbers in this table are for media in ready-to-use packaging.*

Related products

Product	Size	Catalog number
Fibroblast Growth Medium (Ready-to-use)	500 ml	C-23010
Fibroblast Growth Medium Kit	500 ml	C-23110
Fibroblast Basal Medium	500 ml	C-23210
Fibroblast Basal Medium, phenol red-free	500 ml	C-23215
Fibroblast Growth Medium SupplementMix	for 500 ml	C-39315
Fibroblast Growth Medium SupplementPack	for 500 ml	C-39310
Fibroblast Growth Medium 2 (Ready-to-use)	500 ml	C-23020
Fibroblast Growth Medium 2 Kit	500 ml	C-23120
Fibroblast Basal Medium 2	500 ml	C-23220
Fibroblast Basal Medium 2, phenol red-free	500 ml	C-23225
Fibroblast Growth Medium 2 SupplementMix	for 500 ml	C-39325
Fibroblast Growth Medium 2 SupplementPack	for 500 ml	C-39320

Product	Size	Catalog number
Fibroblast Growth Medium 3 (Ready-to-use)	500 ml	C-23025
Fibroblast Growth Medium 3 Kit	500 ml	C-23130
Fibroblast Basal Medium 3	500 ml	C-23230
Fibroblast Basal Medium 3, phenol red-free	500 ml	C-23235
Fibroblast Growth Medium 3 SupplementMix	for 500 ml	C-39345
Fibroblast Growth Medium 3 SupplementPack	for 500 ml	C-39350
DetachKit	30 ml 125 ml 250 ml	C-41200 C-41210 C-41220
Cryo-SFM Plus	30 ml 125 ml	C-29920 C-29922
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
HAoAF Pellet	1 million cells per pellet	C-14037
HCF Pellet	1 million cells per pellet	C-14036
HUF Pellet	1 million cells per pellet	C-14038

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