

Instruction Manual

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

PromoCell offers a range of Human Fibroblasts produced at PromoCell's cell culture facility. The cells are isolated from normal human juvenile foreskin or adult skin from different locations. In addition, PromoCell offers Fibroblasts from adult lung tissue and aortic adventitial tissue.

Shortly after isolation, all PromoCell Human Fibroblasts are cryopreserved at passage 2 (P2) using PromoCell's proprietary, serum-free freezing medium, Cryo-SFM. Each cryo vial contains more

than 500,000 viable cells after thawing. Proliferating cell cultures are made from 500,000 cryopreserved cells that have been thawed and cultured for three days at PromoCell.

Quality Control

Rigid quality control tests are performed for each lot of Human Fibroblasts.

They are tested for cell morphology, adherence rate and cell viability. Flow cytometric analyses of the fibroblast specific antigen CD90 are performed for each lot. Growth performance is tested through multiple passages up to 10 or 15 population doublings (PD) under culture conditions without antibiotics and 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).

A detailed certificate of analysis (CoA) for each lot can be downloaded at: www.promocell.com/coa

Intended Use

PromoCell 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, start immediately with the protocol for cryopreserved cells (see page 2) or the protocol for proliferating cells (see page 3).

Use aseptic techniques and a laminar flow bench.

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.

1. Prepare the medium

Calculate the needed 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 re-tighten. Immerse the vial into a water bath (37°C) just up to 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 prewarmed 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. The cells should be subcultured according to the subcultivation protocol (see page 4) once they have reached 70–90% confluency.



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

Protocol for Proliferating Cells

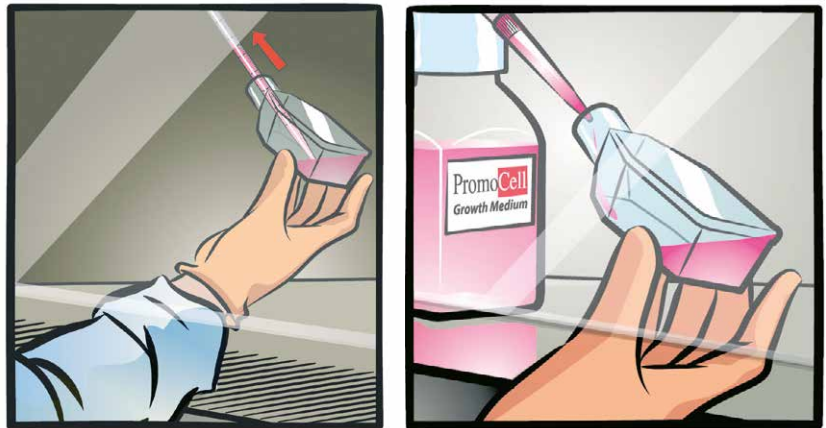
1. Incubate the cells

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 the transportation.



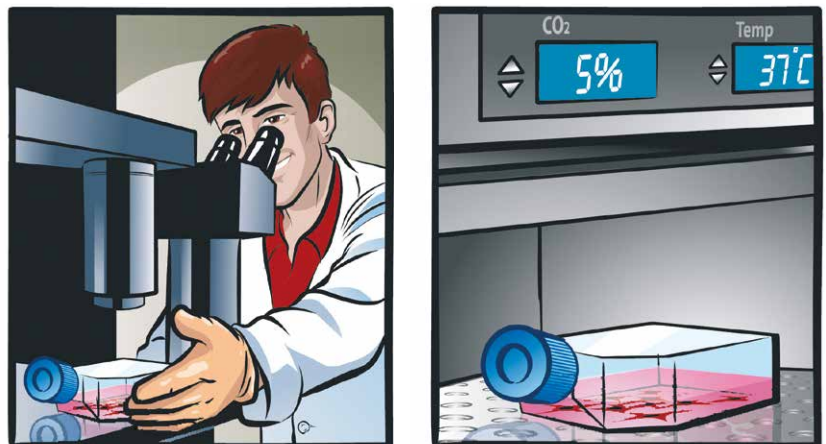
2. Replace the transport medium

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). The cells should be subcultured according to the subcultivation protocol (see page 4) once they have reached 70–90% confluency.

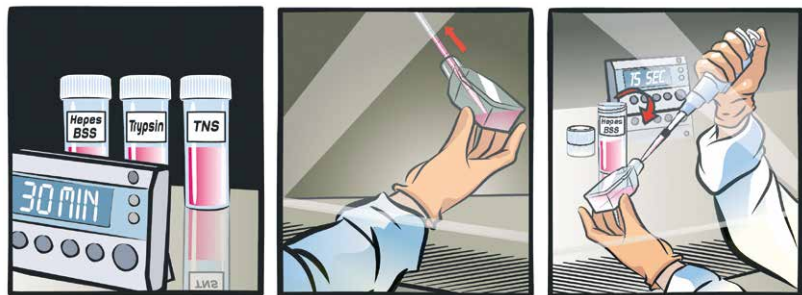


Use aseptic techniques and a laminar flow bench.

Subcultivation Protocol

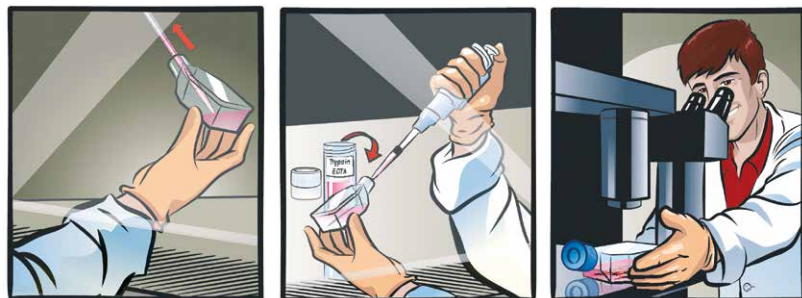
1. Prepare the reagents and wash the cells

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.



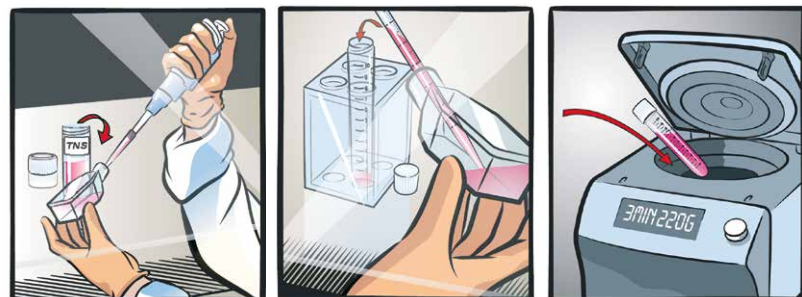
2. Detach the cells

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.



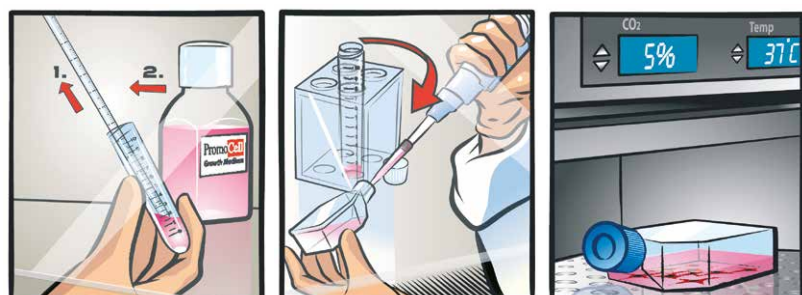
3. Neutralize the trypsin and harvest the cells

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 resuspend 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 Growth Medium. Place the vessels in an incubator (37°C, 5% CO₂) and change the media every two or 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 ⁺ Smooth muscle α -actin ⁺ Slow muscle myosin ⁺	> 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
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	30 ml 125 ml	C-29910 C-29912
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

Email: 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