

Instruction Manual

Product	Size	Catalog Number
Human Osteoblasts (HOB)	500,000 cryopreserved cells	C-12720
	500,000 proliferating cells	C-12760

Product Description

Human osteoblasts are a highly specialized cell type of mesenchymal origin involved in bone formation and remodeling. *In vivo*, they produce the osteoid, an extracellular matrix rich in collagen type I. In the course of their natural maturation process into osteocytes, called osteogenesis, they become embedded in the bone matrix and stop proliferating. PromoCell Human Osteoblasts (HOB) are produced at PromoCell's cell culture facility and are available from knee and hip femoral tissue. Shortly after isolation, all PromoCell Human Osteoblasts 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 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 PromoCell Human Osteoblasts.

They are tested for cell morphology, adherence rate, and cell viability. Furthermore, histochemical tests for alkaline phosphatase and bone mineralization are carried out for each lot (see page 5). Growth performance is tested through multiple passages up to 10 population doublings (PD) under culture conditions lacking antibiotics and antimycotics.

In addition, all cells have been tested for the absence of HIV-1, HIV-2, HBV, HCV, 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 Osteoblasts 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 have to be processed immediately (see page 3).

Start immediately after delivery.

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.



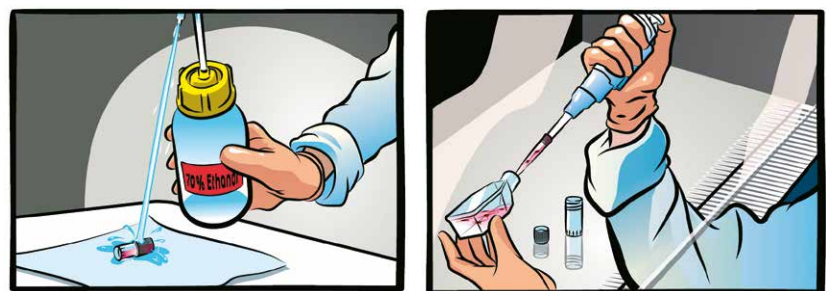
2. Thaw the cells

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 retighten. 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 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.



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

Protocol for Proliferating Cells

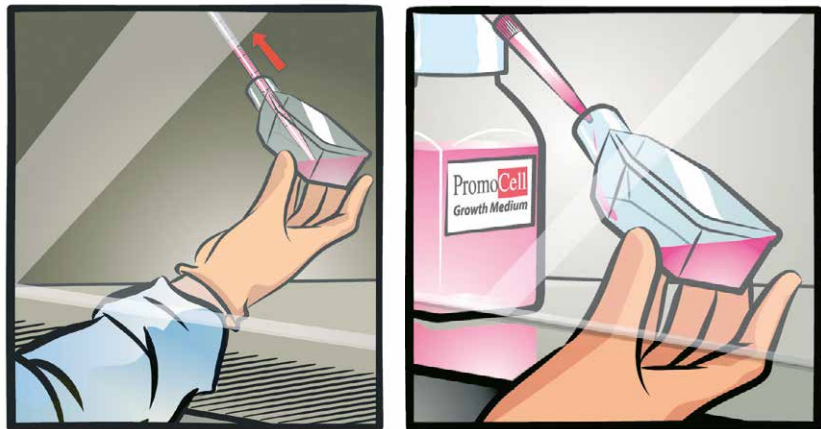
1. Incubate the cells

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.



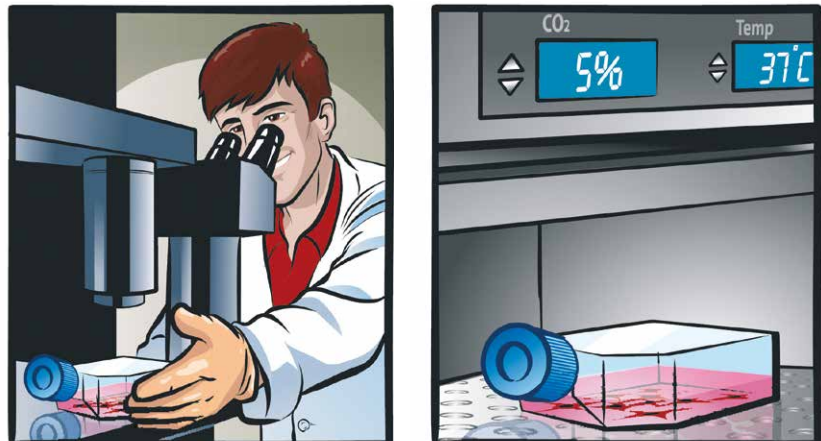
2. Replace the transport medium

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). 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% confluency.

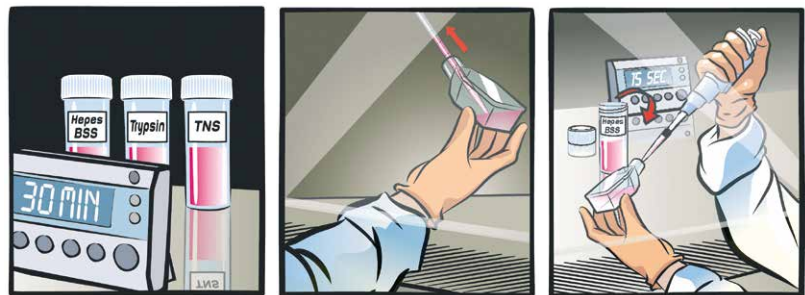


Use aseptic techniques and a laminar flow bench.

Subcultivation Protocol

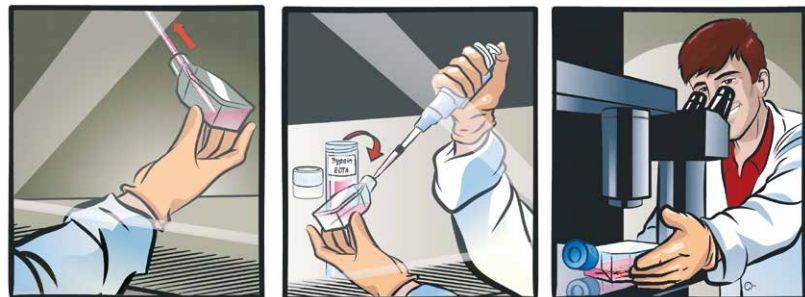
1. Prepare the reagents and wash the cells

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.



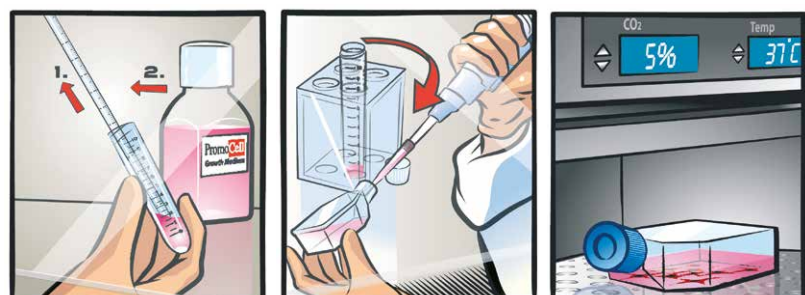
3. Neutralize the trypsin and harvest the cells

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 to three days.



Specifications

Product	Recommended Culture Media*	Plating Density	Passage after Thawing	Marker	Population Doublings
Human Osteoblasts (HOB)	C-27001 C-27020	10,000–20,000 cells per cm ²	P2	Alkaline phosphatase ⁺ Mineralization tested	> 10

Related Products

Product	Size	Catalog Number
Osteoblast Growth Medium (Ready-to-use)	500 ml	C-27001
Osteoblast Basal Medium	500 ml	C-27010
Osteoblast Basal Medium, phenol red-free	500 ml	C-27015
Osteoblast Growth Medium Supplement Mix	for 500 ml	C-39615
Osteoblast Mineralization Medium (Ready-to-use)	100 ml	C-27020
DetachKit	30 ml 125 ml 250 ml	C-41200 C-41210 C-41220
Cryo-SFM	30 ml 125 ml	C-29910 C-29912
HOB Pellet	1 million cells per pellet	C-14071
BMP-2, human, recombinant (HEK)	10 µg	C-67309
sRANKL, human, recombinant (E. coli)	10 µg	C-63200
Vitamin D ELISA Kit, human	10 µg	PK-EL-K2109
TGF-beta 1 ELISA Kit, human	96 tests	PK-EL-63506
Rabbit Anti-Human Osteopontin/SPP-1 (pAB)	100 µg	PK-AB718-7443
PromoFectin cell transfection reagent	0.1 ml 0.5 ml 1 ml	PK-CT-2000-10 PK-CT-2000-50 PK-CT-2000-100
MATra – magnet-assisted cell transfection	Magnet Plate 50 µl MATra-si Reagent	PK-CT-2011-000 PK-CT-2021-005

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

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**