

DetachKit / DetachKit 2

Instruction manual

Product	Size	Catalog number
DetachKit	30 ml	C-41200
	125 ml	C-41210
	250 ml	C-41220
DetachKit 2	30 ml	C-41202
	125 ml	C-41212
	250 ml	C-41222

DetachKit	DetachKit 30 C-41200	DetachKit 125 C-41210	DetachKit 250 C-41220
Contains HEPES BSS Trypsin/EDTA TNS	C-40000 C-41000 C-41100	C-40010 C-41010 C-41110	C-40020 C-41020 C-41120

DetachKit 2	DetachKit 2-30 C-41202	DetachKit 2-125 C-41212	DetachKit 2-250 C-41222
Contains HEPES BSS Trypsin/EDTA 2 TNS	C-40000 C-41002 C-41100	C-40010 C-41012 C-41110	C-40020 C-41022 C-41120

Product description

Our DetachKits were designed for the safe and efficient detachment of primary human cells in routine subculturing. DetachKit 2 is especially valuable for more sensitive cell types.

Each DetachKit consists of three components: HEPES Buffered Saline Solution (BSS), Trypsin/EDTA Solution and Trypsin Neutralizing Solution (TNS).

The DetachKit comes with Trypsin/EDTA (0.04%/0.03%) whereas the DetachKit 2 contains a lower concentration of Trypsin/EDTA (0.025%/0.01%). HEPES BSS contains 30 mM HEPES, D-Glucose, NaCl, KCL and NA-Phosphate. TNS contains 0.05% Trypsin Inhibitor from soybean and 0.1% Bovine Serum Albumin (BSA). All three components

(Trypsin/EDTA, HEPES BSS, and TNS) contain Phenol Red.

Storage and stability

Store at -20°C in the dark after arrival for long term storage. If stored properly, the product is stable until the expiration date stated on the label. The product is shipped with cool packs, so may arrive thawed and can be re-frozen without any loss of quality. However, avoid repeated freeze/thaw cycles. Thawed solutions can be stored at 4 to 8°C and used for up to 6 weeks.

Quality control

All lots of DetachKits are subjected to comprehensive quality control tests. Each lot is

routinely tested for use with primary human cells, the absence of cytotoxicity, and physical parameters such as osmolality and pH level. Approved in-house lots are used as a reference.

In addition, all lots have been tested for the absence of microbial contaminants (fungi, and bacteria).

Intended use

The products are for *in vitro* research use only and not for diagnostic or therapeutic procedures. For safety precautions please see appropriate MSDS.

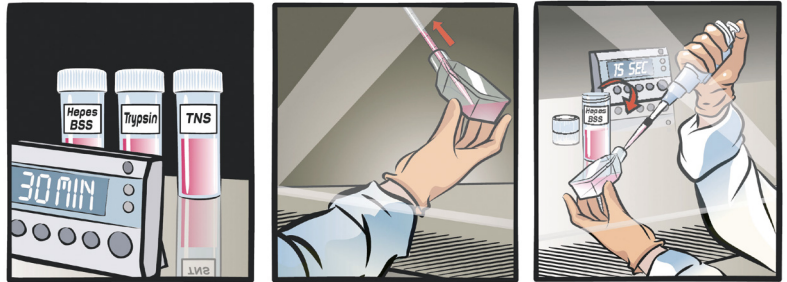
Subcultivation protocol

Use aseptic techniques and a laminar flow bench.

1

Prepare the reagents and wash the cells

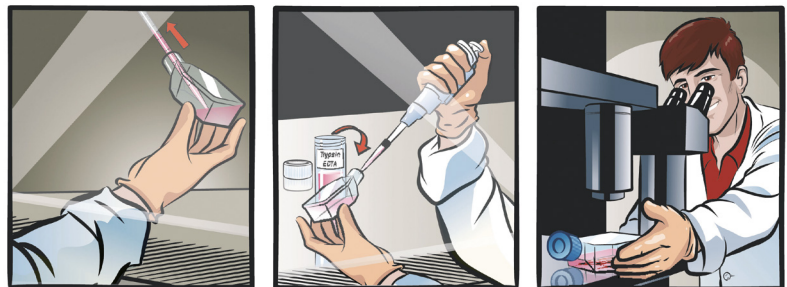
Place the DetachKit at room temperature for at least 30 minutes to adjust the temperature of the reagents. Carefully aspirate the medium from the culture vessel. Add 100 μ l HEPES BSS Solution per cm^2 of vessel surface to wash the cells and agitate the vessel carefully for 15 seconds.



2

Detach the cells

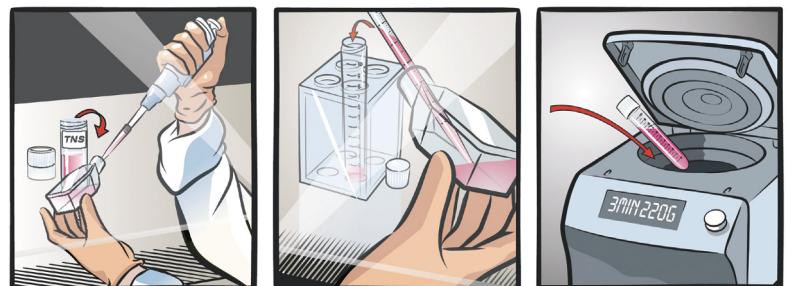
Carefully aspirate the HEPES BSS from the culture vessel. Add 100 μ l Trypsin/EDTA Solution per cm^2 of vessel surface. Note: We recommend detaching the cells at room temperature. Close the vessel and examine the cells under a microscope. When the cells start to detach, gently tap the side of the vessel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



3

Neutralize the trypsin and harvest the cells

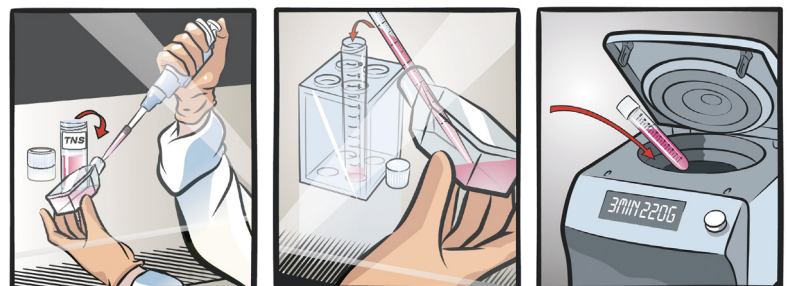
Add 100 μ l Trypsin Neutralizing Solution per cm^2 of vessel surface and gently agitate. Carefully aspirate the cell suspension and transfer it to a centrifugation tube. Spin down the cells for 3 minutes at 220 x g.



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. Determine the cell number and plate the cells according to the recommended seeding density in new cell culture vessels containing PromoCell Cell Growth Medium pre-warmed to 37°C. Place the vessels in an incubator (37°C, 5% CO_2).



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