

Reagents for Cell Detachment

Instruction Manual

Product	Size	Catalog Number
Trypsin/EDTA (0.04% (w/v)/ 0.03% (w/v))	30 ml	C-41000
	125 ml	C-41010
	250 ml	C-41020
Trypsin/EDTA 2 (0.025% (w/v)/ 0.01% (w/v))	30 ml	C-41002
	125 ml	C-41012
	250 ml	C-41022
Trypsin Neutralizing Solution (0.05% (w/v) Trypsin Inhibitor in 0.1% (w/v) BSA)	30 ml	C-41100
	125 ml	C-41110
	250 ml	C-41120

Product Description

Trypsin is still the most widely employed enzyme used for cell detachment. Our Trypsin Solutions exhibit standardized enzyme activity and contain trypsin from extra-pure lots.

Storage and Stability

Store at -20°C in the dark immediately after arrival for long term storage. The product is shipped with cool packs, so may arrive thawed and can be refrozen 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. If stored properly, the product is stable until the expiration date stated on the label.

Quality Control

All lots of PromoCell Cell Detachment Reagents are subjected to comprehensive quality control tests. Each lot is routinely tested for biological function, 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, 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.

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