

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

Description

Trypsin is still the most widely employed enzyme used for cell detachment. PromoCell's 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. Refrigerate thawed solu-

tion at 4 to 8°C. If stored properly, the product is stable until the expiry 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.