

CellCelector Flex

Capillaries, Tips and
Consumables

Simplifying Progress

SARTORIUS

CellCelector Flex Capillaries

A wide range of capillaries and disposable tips are available for the CellCelector Flex for all different application fields. All capillaries and tips are optimized for the specific cell

types and picking modules to guarantee a gentle, precise and contamination-free picking of cells, clusters and colonies.

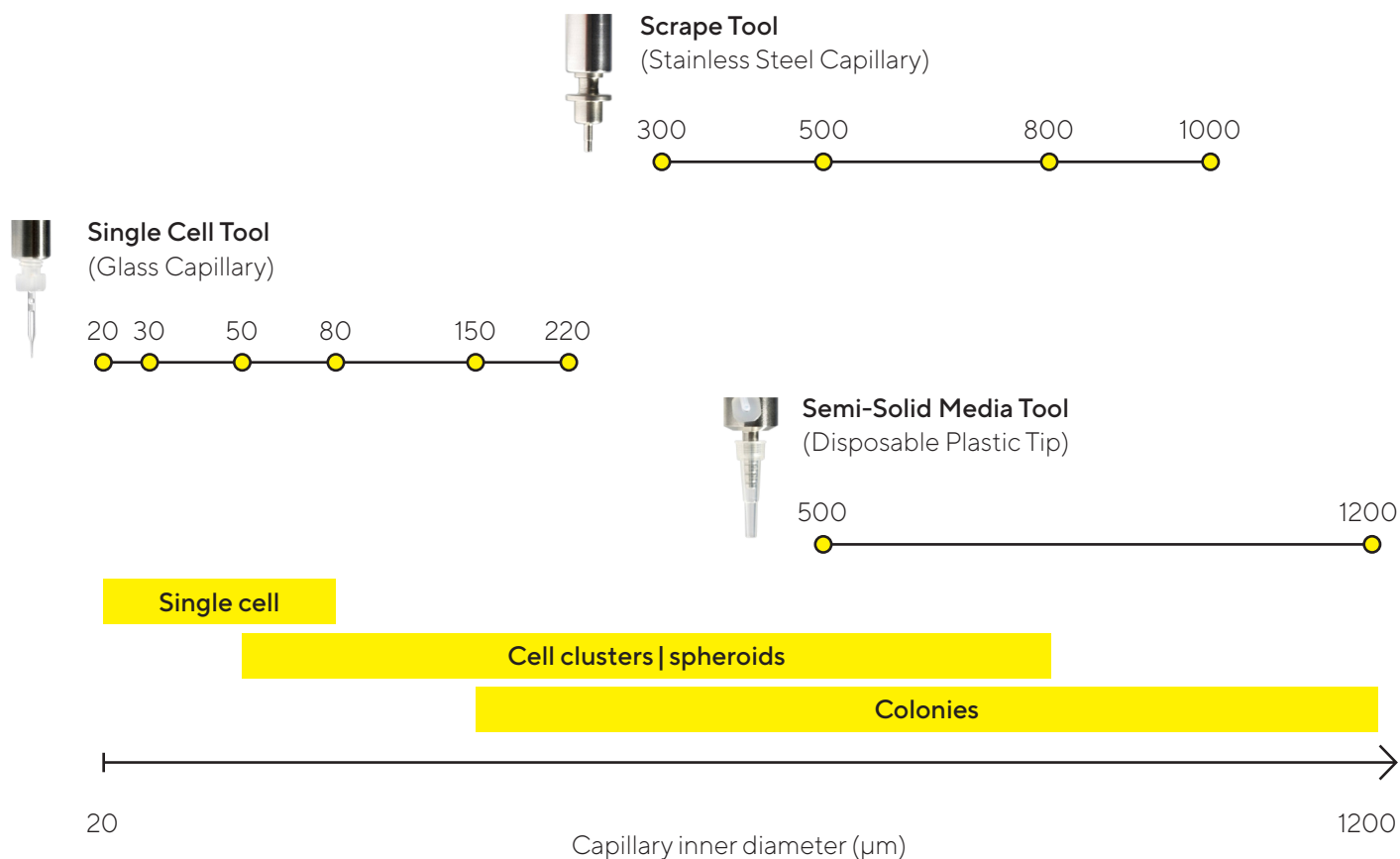


Figure 1

Note. Available tips and capillary diameters for the CellCelector Flex

Disposable Plastic Tips for Semi-Solid Media

The plastic tips are specifically designed for the picking of 3D colonies from semi-solid media like Methylcellulose, Agarose or Matrigel.

The tips are delivered in racks of 96 tips and can be loaded straight onto the CellCelector Flex.



Capillaries for Precise Clone Picking from Semi-solid Media (Picking module CC2003)

Product Number

CellCelector Box PrecisionTip 500, sterilized (96 plastic tips, inner diameter 500 μm)

CC0028

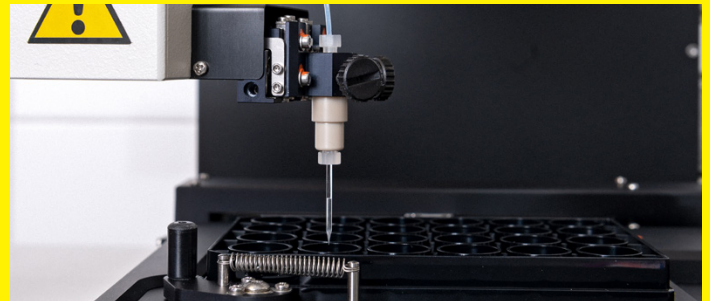
CellCelector Box PrecisionTip 1200, sterilized (96 plastic tips, inner diameter 1200 μm)

CC0027

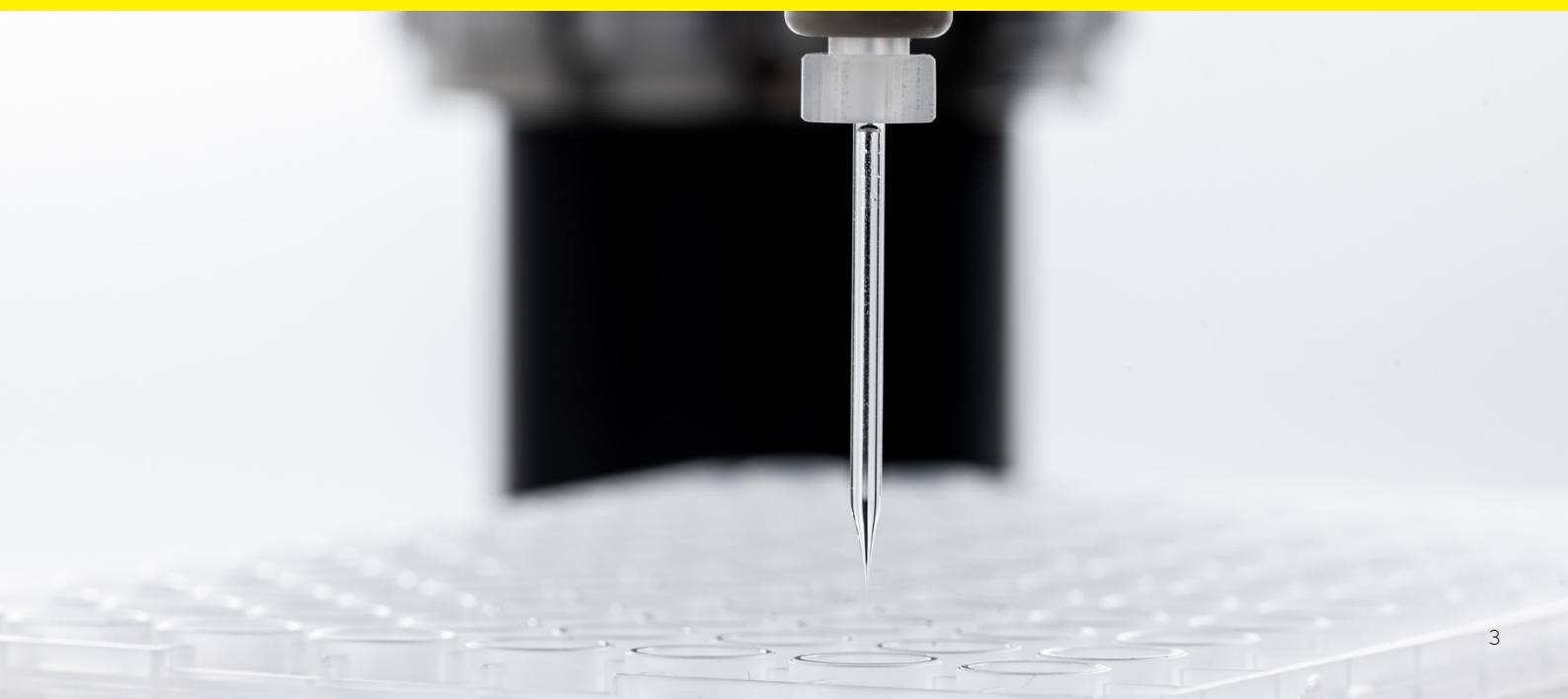
Glass Capillaries for Single Cell Picking and Nanowells

The glass capillaries are designed for the isolation of single cells, clusters, spheroids, organoids and small parts of colonies. Small diameter capillaries allow targeted picking of individual cells without any cross-contamination by adjacent cells.

Larger diameter capillaries enable picking of clusters or spheroids without creating damaging sheer forces keeping the picked cluster intact.

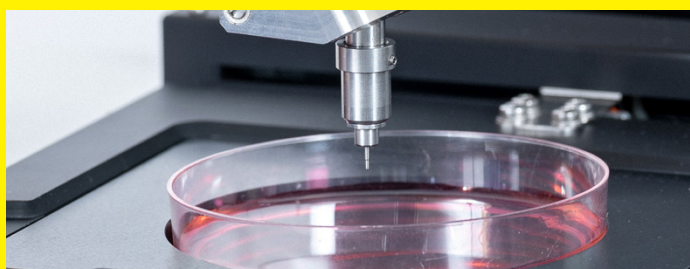


Capillaries for Single Cell Picking (Picking Modules CC2006 and CC2009)	Product Number
CellSelector Single Cell Glass Capillaries, 20 μm (pack of 10, individually packed)	CC0048
CellSelector Single Cell Glass Capillaries, 30 μm (pack of 10, individually packed)	CC0025
CellSelector Single Cell Glass Capillaries, 40 μm (pack of 10, individually packed)	CC0049
CellSelector Single Cell Glass Capillaries, 50 μm (pack of 10, individually packed)	CC0007
CellSelector Single Cell Glass Capillaries, 60 μm (pack of 10, individually packed)	CC0050
CellSelector Single Cell Glass Capillaries, 80 μm (pack of 10, individually packed)	CC0008
CellSelector Single Cell Glass Capillaries, 150 μm (pack of 10, individually packed)	CC0018
CellSelector Single Cell Glass Capillaries, 170 μm (pack of 10, individually packed)	CC0023
CellSelector Single Cell Glass Capillaries, 190 μm (pack of 10, individually packed)	CC0024
CellSelector Single Cell Glass Capillaries, 220 μm (pack of 10, individually packed)	CC0019



Metal Capillaries for Adherent Colonies

The capillaries are designed for the picking of adherent colonies directly from culture vessels without any trypsinization. Made of stainless steel, these capillaries are reusable and autoclavable. Offered in racks of 96 capillaries they can be loaded directly on the deck tray of the CellCelector Flex. Each pack is delivered with a second rack for collection of used capillaries as well as a custom container for autoclaving.



Capillaries for Adherent Colony Picking (Picking Module CC2008)

Product Number

CellCelector Package ScrapeTip 300 (includes a rack of 96 autoclavable metal ScrapeTip, 300 μm , a storage container for autoclaving and a deposit rack)

CC0040

CellCelector Package ScrapeTip 500 (includes a rack of 96 autoclavable metal ScrapeTip, 500 μm , a storage container for autoclaving and a deposit rack)

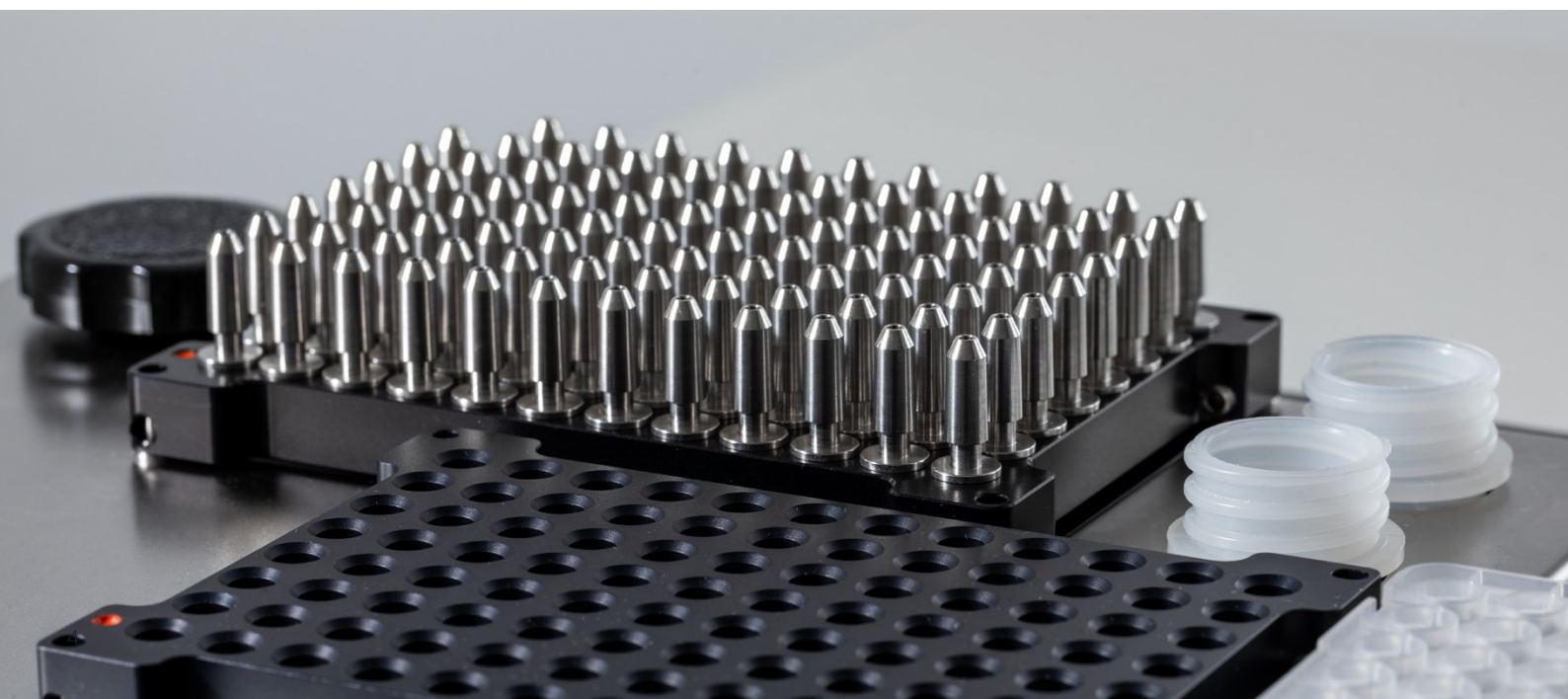
CC0031

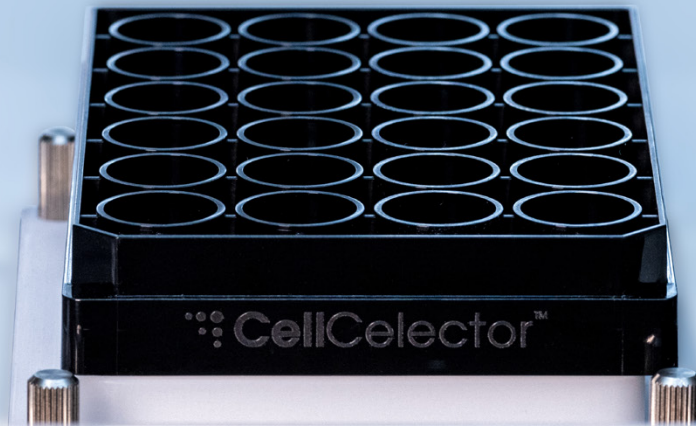
CellCelector Package ScrapeTip 800 (includes a rack of 96 autoclavable metal ScrapeTip, 800 μm , a storage container for autoclaving and a deposit rack)

CC0032

CellCelector Package ScrapeTip 1000 (includes a rack of 96 autoclavable metal ScrapeTip, 1,000 μm , a storage container for autoclaving and a deposit rack)

CC0009





CellCelector Flex Nanowell Arrays

Nanowell arrays for single-cell screening and isolation as well as single cell cloning workflows are easy to use and enable high-throughput and cost-effective alternatives to conventional limiting dilution, FACS and other single cell screening and cloning methods. Our nanowell arrays are

available in various formats from 1 well chamber slides to 6 and 24 well plates. They are featuring from 100,000 to several millions of nanowells per plate | slide at nanowell volumes from several nanoliter down to a few picoliter.

4 nl Nanowell Plates for Single Cell Cloning (S200)

CellCelector Flex S200 nanowell plates are ideal for:

- High-throughput single cell cloning and screening (CHO, HEK, PER.C6 etc.)
- Single cell cloning after gene editing (e.g. stem cells like iPSC etc.)
- Isolation of sets of defined cell numbers (e.g. 1, 5, 20, 50, 100) for reference samples
- Spheroid formation, imaging and screening

Product	Plate S200-100 100K, 24 well, ULA	Plate S200-100 100K, 24 well, Plasma
Product Number	CC0079	CC0075
Qty / Pack	5	5
Number of Wells Plate	24	24
Nanowell Volume	4 nl	4 nl
Nanowell Shape	Square	Square
Nanowell Size	200 µm	200 µm
Nanowell Depth	100 µm	100 µm
Nanowells per Well	3,000	3,000
Nanowells per Plate	72,000	72,000
Nanowell Surface	ULA	Plasma treated
Rec. Glass Capillary Diameter	80 µm	80 µm

Sub-nanoliter Nanowell Arrays for Single Cell Screening and Rare Single Cell Isolation

CellSelector Flex sub-nanoliter nanowell arrays are perfectly suited for:

- Screening of antibody | protein producing cells (e.g. B-Cells) through secretion assays
- Co-location of different cell types in the same nanowells for functional studies or receptor | reporter cell assays
- Isolation and analysis of rare single cells like circulating tumor cells and fetal cells
- Hybridoma single cell cloning and screening

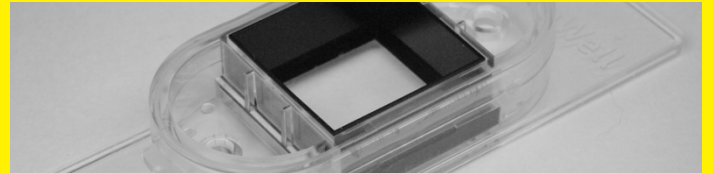
Product	Nanowell plate H100-100 330K, 24 well	Nanowell plate H100-100 360K, 6 well	Nanowell plate U40-40 850K, 24 well	Nanowell plate U40-40 1M, 6 well	Nanowell slide U40-40 200K, 2 well	Nanowell slide U25-25 400K, 2 well
Product Number	CC0084	CC0080	CC0096	CC0097	CC0083	CC0081
Qty Pack	5	5	5	5	5	5
Number of Wells/Plate	24	6	24	6	2	2
Nanowell Volume	0.9 nl	0.9 nl	50 pl	50 pl	50 pl	12 pl
Nanowell Shape	Hex	Hex	UFO	UFO	UFO	UFO
Nanowell Size	100 µm	100 µm	40 µm	40 µm	40 µm	25 µm
Nanowell Depth	100 µm	100 µm	40 µm	40 µm	40 µm	25 µm
Nanowells per Well	14,000	60,000	36,000	160,000	50,000	200,000
Nanowells per Plate	330,000	360,000	850,000	1,000,000	200,000	400,000
Nanowell Surface	Non-treated	Non-treated	Non-treated	Non-treated	Non-treated	Non-treated
Rec. Glass Capillary Diameter	50 µm	50 µm	30 µm	30 µm	30 µm	20 µm



SIEVEWELL Nanowell Arrays for Rare Single Cell Isolation Workflows

The SIEVEWELL chip consists of a thin membrane with nanowells. The outer dimensions of the chip correspond to the standard microscope slide format. Within each nanowell there are 2 pores with a diameter of 2 μm . The hexagonal shape of the wells is ideal for automated cell detection and cell counting. Due to its unique sieve structure SIEVEWELL arrays are a simple and perfect solution for:

- On-chip staining of rare single cells, like circulating tumor cells, fetal cells etc. without any cell loss
- Single cell capture from cell suspensions: just pipetting without pumps.



Product	SIEVEWELL Slide 370K	SIEVEWELL Slide 90K
Product Number	CC0085	CC0098
Qty Pack	5	5
Number of Wells Slide	1	1
Nanowell Volume	8.6 μl	108 μl
Nanowell Shape	Hex	Hex
Nanowell Size	20 μm	50 μm
Nanowell Depth	25 μm	50 μm
Nanowells per Slide	370,000	90,000
Nanowell Surface	Non-treated	Non-treated
Rec. Glass Capillary Diameter	20 μm	20 μm

CellCelector Flex MagnetPick Slides

The MagnetPick slides are designed for the specific isolation and picking of rare single cells (like CTCs or fetal cells) from suspension that have been processed utilizing positive immuno-magnetic enrichment methods (e.g. CellSearch[®], Isoflux[™], MagSweeper[™]).

They can only be used in combination with the MagnetPick adapter.


Description	Product Number
MagnetPick Slides 300 1 pack of 72 MagnetPick glass slides for MagnetPick 300 adapter; Each slide contains one hydrophilic sample area for sample volumes of up to 300 μl	CC0059

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For additional information,
visit www.sartorius.com