

## Drug Testing & Cell and Gene Therapy

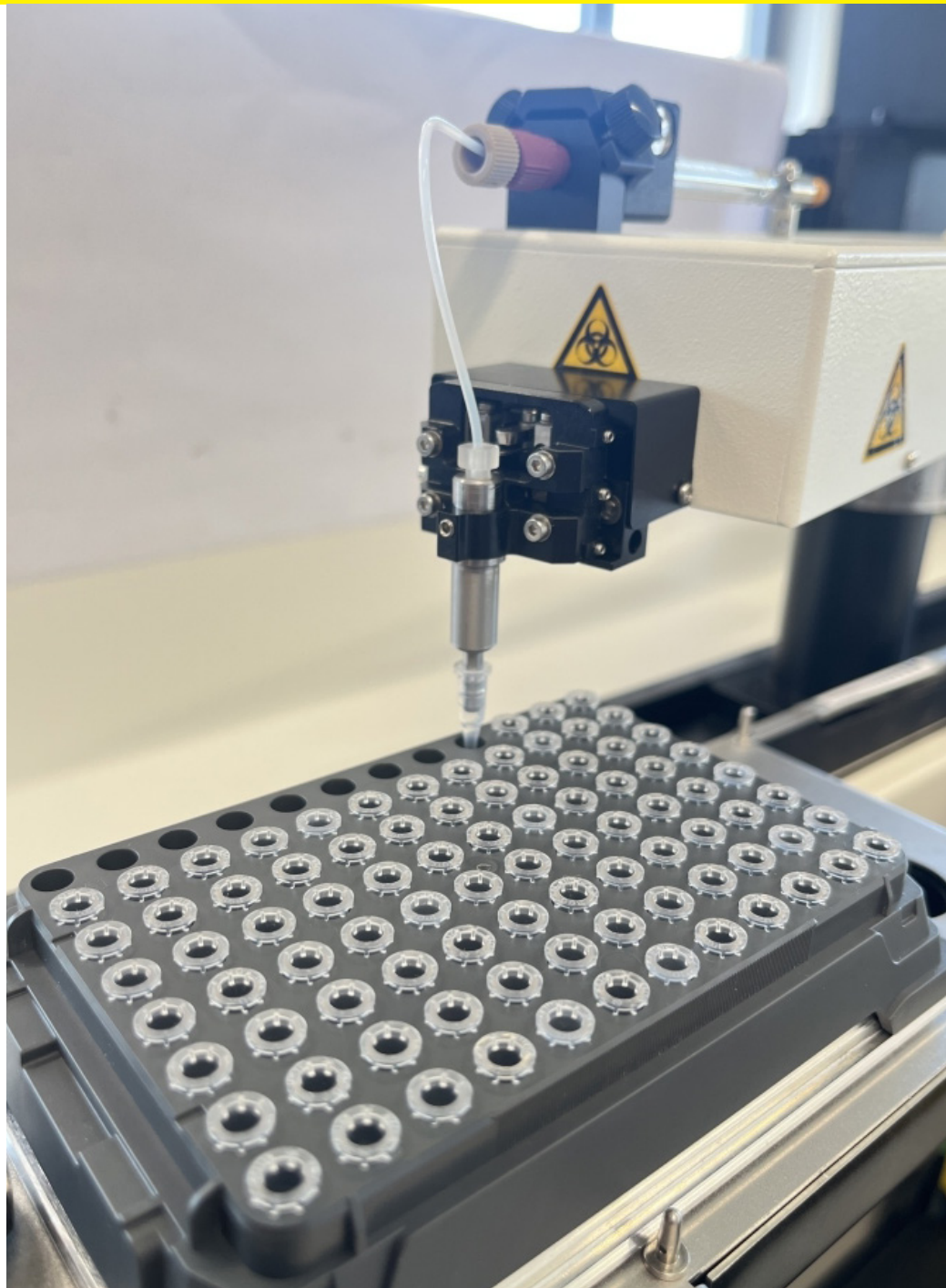
High Performance Precision Tips for the CellCelector Flex

### The Ideal Combination for Your Cell Selection and Retrieval

- Highly accurate and gentle cell isolation and transfer, with excellent post-picking viability rates
- Automated and manual selection of clones, colonies and objects of interest
- Full documentation of the picking process, including both pre- and post-picking image verification
- Contamination-free automated picking
- Validated liquid and residual-free tip deposition to ensure full content transfer

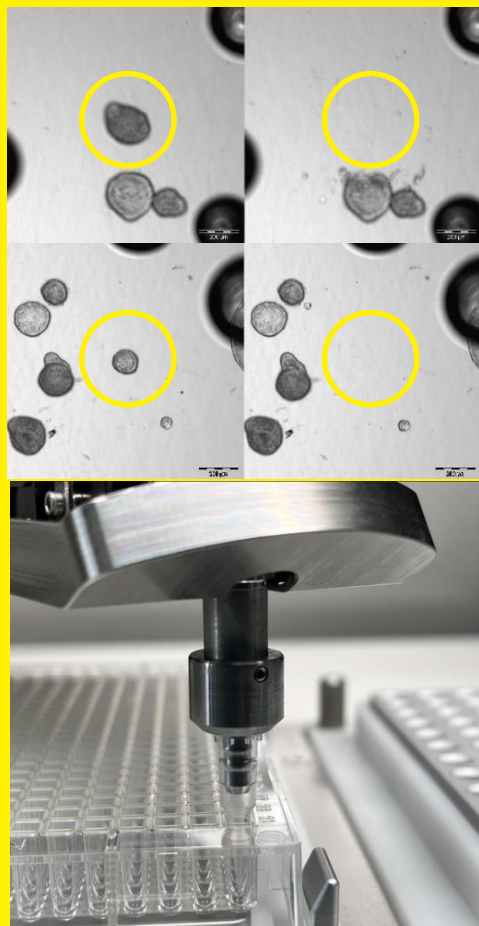


Contact us for efficient  
cell selection and retrieval:  
[CellCelector\\_Support@Sartorius.com](mailto:CellCelector_Support@Sartorius.com)



# High Performance Precision Tip Features

High Performance Precision (HPP) Tips were developed to facilitate the automatic exchange of disposable plastic tips to ensure contamination-free cell selection and transfer. Combined with the CellCelector Flex, HPP Tips provide researchers with a gentle and accurate transfer of clones, organoids, spheroids and adherent biological objects for all drug discovery and cell and gene therapy applications, resulting in continually high post-picking viability rates to help facilitate further analysis.

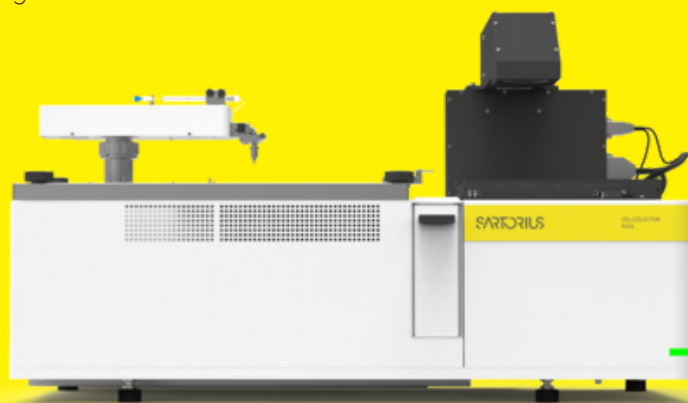


## Picking Advantages

- Flexible range of tip sizes, including 300  $\mu\text{m}$ , 800  $\mu\text{m}$  and 1500  $\mu\text{m}$  diameters
- Fully automated HPP Tip pick-up and release
- Accurate and reproducible pick-up positions following multiple object transfer and tip exchange
- Increased available area for cell picking compared to the stainless steel scrape tips, including 6 well plates (83% vs. 73% of well surface), 24 well plates (63% vs. 40% of well surface) and Petri dishes (89% vs. 82% of well surface)
- Validated organoid picking from a range of Matrigel™ concentrations of up to 100%

## Deposition Advantages

- Validated use of a range of deposition plates, from 6 well plates to 384 well plates
- No liquid residues remaining within HPP Tips, to ensure that all tip contents following aspiration are accurately deposited
- Validated cell viability of up to 100% after 3 days following cell transfer



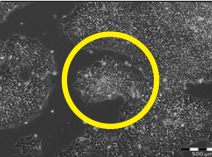
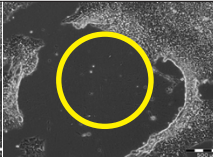
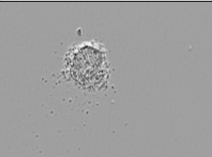
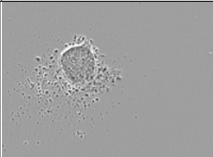
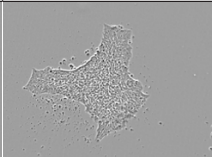
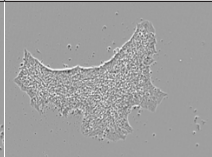
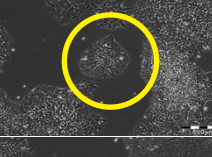
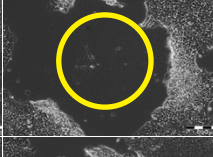
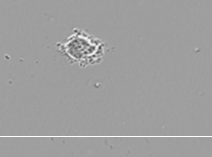
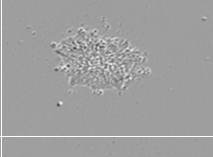
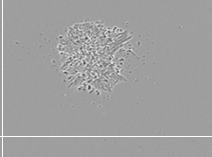
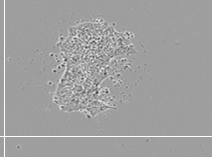

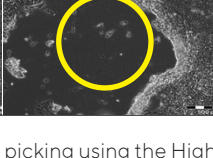
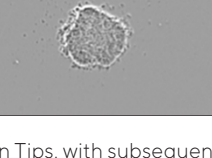


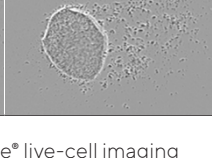
## Ordering Information

Product Number	Description	Pack Size	
CC0002	300 $\mu\text{m}$ High Performance Precision Tips	1 box of 96 HPP Tips	
CC0003	500 $\mu\text{m}$ High Performance Precision Tips	1 box of 96 HPP Tips	Coming Soon
CC0004	800 $\mu\text{m}$ High Performance Precision Tips	1 box of 96 HPP Tips	
CC0005	1000 $\mu\text{m}$ High Performance Precision Tips	1 box of 96 HPP Tips	Coming Soon
CC0006	1500 $\mu\text{m}$ High Performance Precision Tips	1 box of 96 HPP Tips	
CC0011	300 $\mu\text{m}$ High Performance Precision Tips	Pack of 5 boxes of 96 HPP Tips each	Coming Soon
CC0012	500 $\mu\text{m}$ High Performance Precision Tips	Pack of 5 boxes of 96 HPP Tips each	Coming Soon
CC0013	800 $\mu\text{m}$ High Performance Precision Tips	Pack of 5 boxes of 96 HPP Tips each	Coming Soon
CC0014	1000 $\mu\text{m}$ High Performance Precision Tips	Pack of 5 boxes of 96 HPP Tips each	Coming Soon
CC0015	1500 $\mu\text{m}$ High Performance Precision Tips	Pack of 5 boxes of 96 HPP Tips each	Coming Soon
CC2007	Next Generation Precision Tip Picking Module	1 Picking Arm for Scrape and Semi-Solid Media Applications	
CC3010	HPP Tip Deck Tray Adapter	1 Adapter	

# High Performance Precision Tips : Example Results

Adherent induced Pluripotent Stem Cells (iPSCs) were seeded into two wells of a 6-well plate on a CellCelector Flex, before being scanned using a 10x objective in phase contrast. A 96-well cell culture target plate was prepared and the appropriate picking parameters selected. During automated cell transfer, the system recorded before and after sample images to ensure full documentation of the picking process. To further investigate the growth characteristics of the picked iPSCs, the transferred cells were subsequently cultured and imaged using an Incucyte® live-cell analysis instrument for a further 3 days.

A defined volume of medium was aspirated from the target 96-well plate, before iPSC colonies were identified, scraped, and deposited back into designated wells of the target plate. Results highlight a picking efficiency of 100% and a transfer efficiency of 100% for cellular deposition. IPSC growth and associated colony viability was monitored by scanning the plates at two-hour intervals for a period of three days using the Incucyte®, which found a viable colony efficiency of 94%.

CellCelector Images		Destination Well	Incucyte® Image Verification (4x scanning)			
Before Picking	After Picking		Day 0	Day 1	Day 2	Day 3
		C01				
		C02				
		C03				

Adherent iPSC colony picking using the High Performance Precision Tips, with subsequent viable colony assessment using the Incucyte® live-cell imaging platform.

## Germany

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

## USA

Sartorius Corporation  
3874 Research Park Dr.  
Ann Arbor, MI 48108  
Phone +1 734 769 1600



For further information, visit  
[sartorius.com](https://www.sartorius.com)