

# CellCelector

FULLY AUTOMATED CELL PICKING SYSTEM

## HT-NIC: HIGH-THROUGHPUT NANOWELL-BASED IMAGE-VERIFIED CLONING METHOD



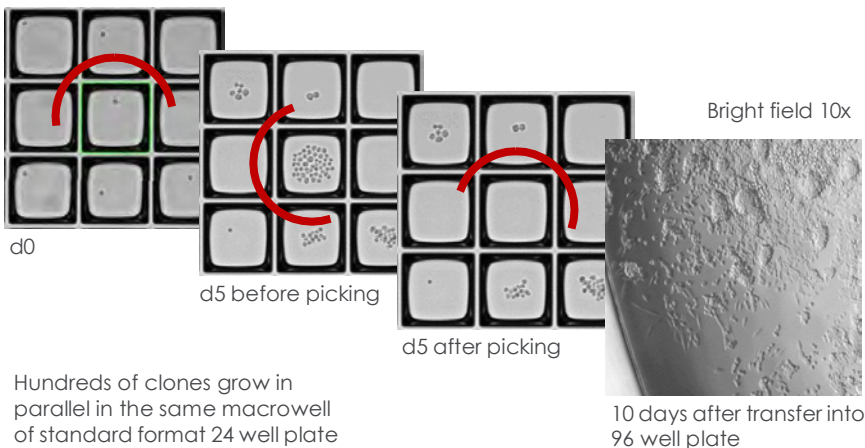
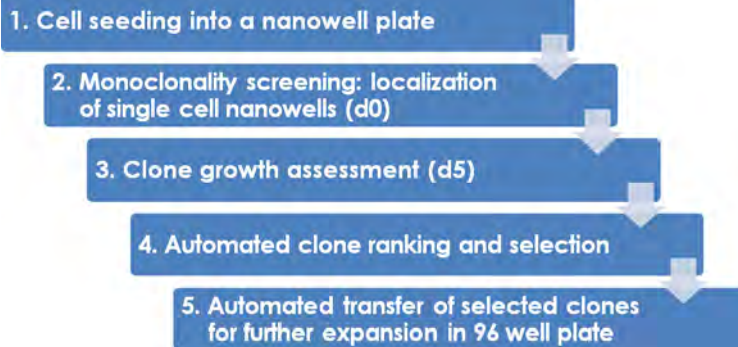
**fast** cell line development  
integrated **monoclonality proof**  
important **time and cost savings**

The ALS CellCelector single cell and colony picking platform can now be used for high-throughput single cell cloning allowing fast generation of clonal production cell lines with one cloning round while providing in-process image-verified monoclonality proof. With integrated monoclonality and clone viability assessment as well as high outgrowth rate after clone transfer to 96 well plates, the CellCelector HT-NIC technology represents an advantageous alternative to limiting dilution or FACS single cell sorting techniques. The method has been developed in collaboration with ProBioGen AG.

  
Automated • Lab • Solutions  
Now part of Sartorius

## Nanowell-Based Single Cell Cloning Workflow

The method uses standard-format 24 well CellCelector Nanowell plates (available from ALS) with thousands of tiny nanowells on the bottom of each well. Bright field scanning of seeded wells followed by the automated identification of all single cell nanowells provides a robust and documented image-based mono-clonality proof. After such mono-clonality scan cells are let to grow several days in an incubator. As early as on Day 4 after seeding the nanowell plate is scanned again and the viable clones resulting from single cells are automatically selected and transferred into 96 or 384 well plates for further growth and productivity assessment.



## Benefits and Key Features

- Faster CLD times (by 5 to 9 weeks)
- Integrated image-verified mono-clonality proof
- High image quality for robust automated label-free single cell detection
- Selection of clones by outgrowth and/or by fluorescence
- 100% selective clone recovery without cross contamination
- High outgrowth efficiency of transferred clones
- Significant cost savings on consumables, media and incubator storage space. **Just one plate per cloning experiment**
- Easy-to-change disposable single-use capillaries
- No routine maintenance necessary

### For More Information

ALS Automated Lab Solutions, now part of Sartorius  
Otto Eppenstein-Str. 30, 07745 Jena, Germany  
Phone: +49 (0) 3641 4820-0  
Fax: +49 (0) 3641 4820-11  
Email: info@als-jena.com

### Germany

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

### USA

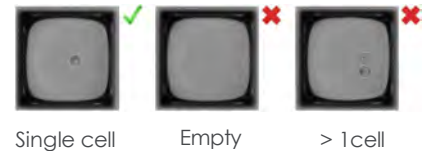
Sartorius Corporation  
565 Johnson Avenue  
Bohemia, NY 11716  
Phone +1 631 254 4249  
Toll-free +1 800 635 2906



ALS CellCelector installed in ALS FlowBox Incubator with temperature, humidity, and CO<sub>2</sub> control. The system can also be placed within a standard biosafety cabinet.

## Mono-clonality Proof

Cells are clearly visible within 4 nL nanowells and can be reliably detected.



## Other Applications

ALS CellCelector is an open platform which can be used for multiple applications:

- Single cell isolation
- scRNA-Seq
- CRISPR/CAS9
- Rare cell detection
- Stem cells
- Semi-solid media cloning (additional module)