

SARTORIUS



Detection, Selection and Isolation of Single Cells and Adherent Colonies

CellCelector™ Flex

Versatile Automated Single Cell and Colony Picking System

The world's most flexible cell picking system

- Unique and fully automated cell imaging and picking system, combining an automated inverted fluorescence microscope, a high-speed scanning stage, and a high-precision microfluidic cell picking robot
- Developed for the detection, selection and isolation of single cells, clusters, spheroids, organoids, single cell clones and adherent colonies
- Used across a variety of research areas, such as circulating tumor cell (CTC) screening, stem cell research, cell line development (CLD) and antibody discovery

Automated Lab Solutions: now part of Sartorius

A Versatile Automated Cell and Colony Picking System

The CellCelector™ single cell and colony picking platform is based on a multi-patented, unique cell and colony isolation technology which provides unmatched flexibility, high-quality results, operational reliability, cost-effectiveness and simple use.

The cell picking technique is predicated on a non-invasive mechanical aspiration process where cells of interest can be detected, identified and selected in a fully automated manner, using either label-free size and morphology characteristics, or fluorescence markers.

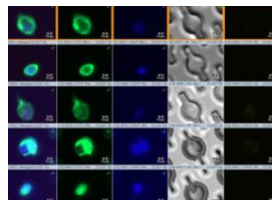
The entire workflow is fully documented by providing live images before, during and after cell picking, as well as cell tracking data from source to destination to ensure full traceability of the screening and cell recovery process.

Key features and specifications of the CellCelector™ include:

- Open platform allowing the development of your own imaging, detection and picking protocols
- Versatility: single cell picking, isolation of clusters, spheroids, organoids, adherent colonies, and single cell clones from a variety of liquid and semi-solid medias
- Up to 100% picking/transfer efficiency
- Extremely gentle cell transfer resulting in high cell integrity and outgrowth rates after cell transfer
- Isolation of living or fixed cells

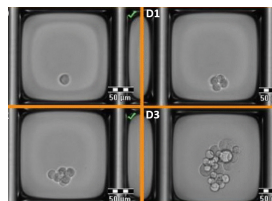
Validated Applications

Single Cell Isolation



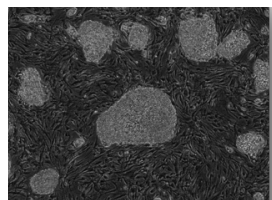
- CTC isolation and analysis
- Rare single cells
- Fetal cells cbNIPT
- Sperm cell isolation
- Protoplasts plant cells
- Single cell heterogeneity
- CRISPR single cell cloning

Bioprocessing



- Hybridoma and CHO clone picking
- Single cell cloning for CLD
- Automatic transfer of bacterial colonies
- 1-day single plasma B cell secretion assays

Stem Cells



- Clonal stem cell picking
- Hematopoietic stem cell colonies
- HSC daughter cell splitting

Learn More:



www.als-jena.com