



Automated • Lab • Solutions
Now part of Sartorius

Time for
Science



 **CellCollector**

Laminar Flow Cabinet for

Working with primary cells and tissue

Flexible and long-term analysis of cells and colonies

Experiments under physiological and sterile conditions



Technical Specifications



Stable and light frame

- WxDxH 1.10m x 0.85m x 2.30m
- Weight: 175 kg / 385 lb
- CellCelector is placed inside
- Height with floor cupboard 2.30 m / 7.5 ft
- Rolls allow for flexible placing in the lab

Filter system

- Filter HEPA H14 (0.1 – 0.3 µm)
- A constant air stream is surrounding the CellCelector with a stream of 0.5 m/sec
- Outgoing air is filtered to avoid room contamination

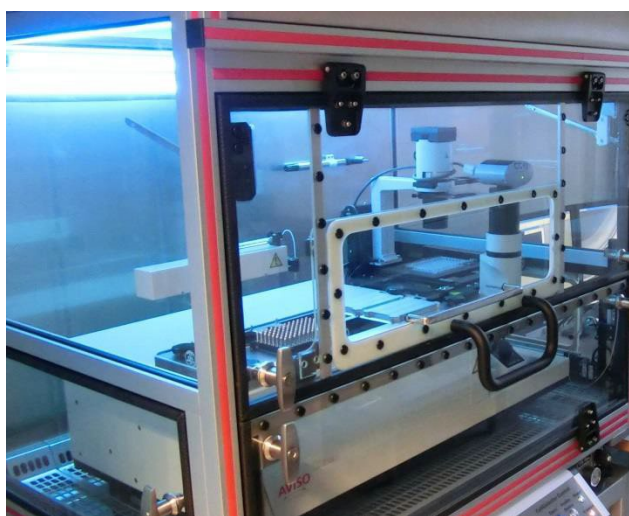


Adjustable temperature up to 40°C

- Individual setting in the control panel
- Temperature control permanently regulates to maintain the set temperature

High CO₂ atmosphere (5%)

- The CO₂ gas flow is permanently regulated and holds constantly at the set level



High humidity atmosphere up to 75%

- The humidity level can be set by the operator on the control panel

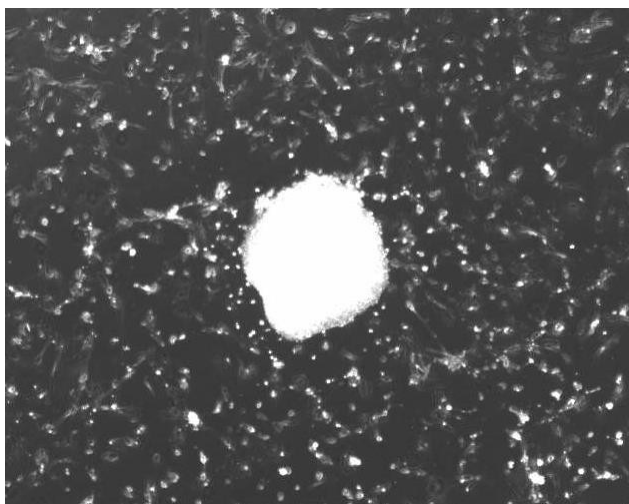
UV-C Illumination

- Provides sterile surfaces
- Four high energy UV-C lamps 4 x 15

Security management system

- Access doors on each side
- Security management system stops all mechanical activity inside the Flowbox once user intervention is necessary

Physiological Conditions for Cell Handling



High CO₂-atmosphere and constant high temperature (37°C) allows for cell selection, analysis, and maintenance of cell lines with the CellCelector.

Advanced level of security against contamination of valuable cell culture material with airborne pathogens. Experiments without time limitations and less differentiation or cell death due to temperature changes and high air Oxygen (hypoxia stress).



Investigation of primary cells not adapted to air Oxygen content (e.g. monocytes, cancer stem cells) and tissue under physiological conditions can reveal new data and models. Remote handling of cells with the CellCelector inside the Flowbox reduces the human factor and sets standards for experimental procedures.



Additional technical equipment and power supply for the CellCelector can easily be stored underneath the Flowbox in the floor cupboard. It is mounted on rolls which allows for flexible placement in the lab room.



Automated • Lab • Solutions
Now part of Sartorius

Time ^{for} Science



ALS Automated Lab Solutions headquarters in Jena, Germany

ALS Automated Lab Solutions GmbH is located in Jena, a dynamic city famous for microscopy and material science. ALS is a specialist for the development of innovative technological solutions for cell biology research and molecular biology. We lift cell culture to a new level of choice and control on the leading edge in cell biology, cell therapy research, regenerative medicine and drug discovery. With automation and standardization of laborious manual procedures, ALS supports science and research for more efficiency and the creation of new methods for the science of tomorrow.

SARTORIUS

For More Information

Jens Eberhardt
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

For additional information,
visit www.sartorius.com

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