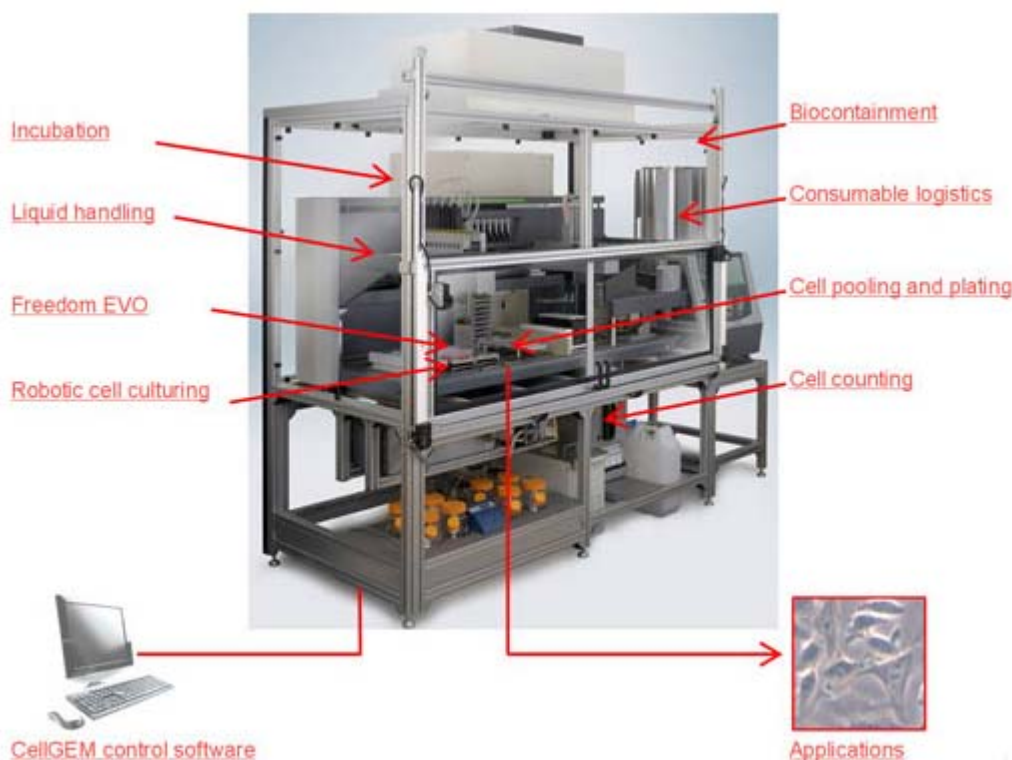


Cellerity™

Consistent production and delivery of ready-to-use cells

Reference 10001BO100398



Fully automated cell culture of adherent cell lines, cell counting, culture maintenance, process scheduling software and hardware, consumable logistics

Cellerity is a modular system which allows standard cell biology processes to be performed for adherent cell lines, including cell feeding, harvesting and passaging in microplate-sized, automation-friendly cell culturing flasks. Tecan suggests using Corning® RoboFlasks™. The Cellerity system can also be used to prepare cells for downstream applications such as cell-based screening or protein expression studies and allows processing of several cell lines in parallel on the system.

CellGEM™ (Cell Growth Expansion and Maintenance) software is a part of the Cellerity system. It is an intuitive software interface that guides users through the process of initiating a cell line expansion request. The CellGEM scheduler calculates optimal times for cell maintenance and production requests. CellGEM automatically instructs the instrument to perform the required actions by using Tecan's Freedom EVOware® Plus pipetting software. CellGEM also maintains a record of the cells grown on the system, including growth rates and passage number. Additionally, CellGEM keeps track of current volumes of media and consumables available in the system and warns users of depleting resources.

The modular Cellerity system can be expanded with a centrifuge allowing suspension cell culturing, or with detection devices to control cell culture confluency.

Scientific instrumentation; not for use in human, clinical or diagnostic procedures. All systems or products are tailor-made one-offs that were exclusively developed on customer demand. All displayed pictures and images originate from development, test or installation phases and do not claim to represent the final installation at customer site.

Tecan and Freedom EVOware are in major countries registered trademarks, and Cellerity and CellGEM are trademarks of Tecan Group Ltd., Männedorf, Switzerland. Corning and RoboFlask are registered trademarks of Corning Incorporated, Corning, USA.
© 2009, Tecan Trading AG, Switzerland, all rights reserved.