

Cell Biology

Monoclonal Antibody Production, Cell Expansion

This system is used for the automated cell expansion and confluence measurement of cell lines which produce monoclonal antibodies. Part of the Monash system.

Technical Reference 30032250



Input

Cells in 96 well microplates

- Cells are provided in 96 well plates

Process

Expansion of cells from 96 well to 24 well and to 6 well plates

- A hit picking application picks the cells marked as hits from the wells of the 96 well plates into the 24 well plates
- The cells are then incubated in the Cytomat@ 48C
- Confluence of the cells in the 24 well plates is measured
- If the threshold is reached the cells are expanded from the 24 well plates to 6 well plates (hit picking application)
- If the threshold is not reached the cells are grown for another day and confluence measurement is repeated
- Again, cells that have reached the growth threshold are expanded to 6 well plates

Output

Cells in 24 well plates

- Cells in 24 well plates, one cell line per plate

Cells in 6 well plates

- Cells in 6 well plates, one cell line per plate

Throughput

- Either 24 x 24 well plates or 96 x 6 well plates per working day

Highlights

Integration of the CloneSelect Imager (Provided by Genetix)

The CloneSelect Imager is used for confluence measurement in 24 well plates and is placed onto an adapter plate in the right front angle of the Freedom EVO®.



Integration of Cytomat 48

The transfer rail of the Cytomat presents the plate directly on the Freedom EVO worktable. Pipetting into the plate can be performed directly on the transfer slide (no ROMA movement necessary).



Scientific instrumentation; not for use in human, clinical or diagnostic procedure. 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, Freedom EVO are in major countries registered trademarks and MCA is a trademark of Tecan Group Ltd., Männedorf, Switzerland. IKA is a registered trademark of IKA Werke GmbH, Hilden, Germany. Cytomat is a trademark of Kendro Laboratory Products GmbH, Hanau, Germany.
© 2008, Tecan Trading AG, Switzerland, all rights reserved.