

Freedom EVOware®

Simple control of complex sample pipetting procedures

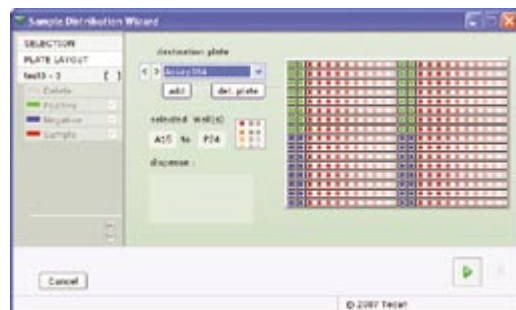
The latest Freedom EVOware® Sample Oriented software, version 1.2, is now available for the Freedom EVO® family of workstations. This software enhances Freedom EVOware Standard and Freedom EVOware Plus with a number of sample-based pipetting, diluting and pooling functions. Freedom EVOware Sample Oriented is ideal for laboratories that need to perform complex pipetting and dilution sequences, and for those with variable joblists, assays and plate definitions. Freedom EVOware Sample Oriented is essential for pooling samples* with the Freedom EVO Clinical and Te-PoolSafe™, and controls primary pooling, sub-pooling, cross pooling and pipetting to archive plates.

The plug-in wizards provide simple, graphically-guided operation for many tasks, including setting up assay plates, and defining pipetting sequences and pre- or serial dilutions. The wizards also manage sample replicates, control reagent addition and volume tracking,

and allow parallel pipetting of samples to one or more assays. The optional loading interface detects whether samples have been manipulated prior to pipetting and assists the user with loading of consumables onto the workstation.

Positive identification of samples, plates and liquids can be carried out using PosID™ 2 or PosID 3 on the Freedom EVO 100, 150 and 200, or using a hand-held scanner on the Freedom EVO 75; sample identities can be tracked throughout pipetting steps and output files created. Freedom EVOware Sample Tracking is available as an optional add-on that creates printed reports for each plate (Sample Tracking Embedded) and provides long-term data storage (Sample Tracking Standard).

Freedom EVOware Sample Oriented can be used with Magellan™, Tecan's sophisticated data processing software for detection instruments, and has LIMS and ASTM E1381/E1392 connectivity.



The software package is included with the Freedom EVO 75, and can be ordered in addition to Freedom EVOware Standard or Plus for other Freedom EVO platforms. Freedom EVOware Sample Oriented allows a 30-day evaluation upon installation of Freedom EVOware on an instrument, before a dongle is required for continued operation.

** Freedom EVO Clinical, pooling and Te-PoolSafe option not available in the USA*

For more information on Tecan's Freedom EVOware Sample Oriented, visit www.tecan.com/freedomevowaresampleoriented

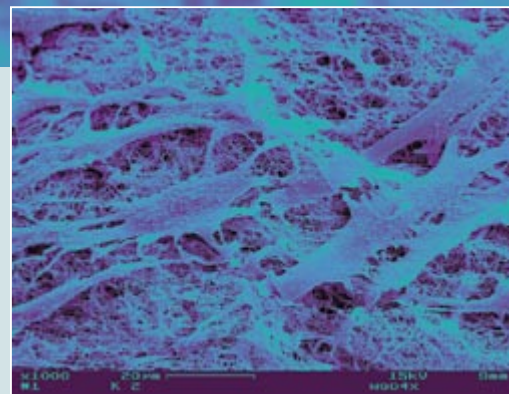
Making advances in regenerative medicine

Tecan is collaborating with tissue engineering researchers at the University of Zurich to develop a novel, automated system for propagation of quality controlled cells that will be suitable for use in regenerative medicine therapies, such as intervertebral disk implants. The project is funded by the CTI, the Swiss Confederation's innovation promotion agency, which supports the transfer of knowledge and technology between universities and businesses.

The new solution will control all necessary cell culture steps, including the isolation,

seeding, proliferation, harvesting and analysis of cell lines and primary cells. Quality control will be maintained through two integrated detection devices that will determine the degree of cell confluence and detect the expression of cell type-specific genetic markers. The platform will be easily adjusted for culture of multiple cell types, providing standardized and affordable treatment procedures for regenerative medicine.

Research application, not for clinical diagnostic use.



SEM (scanning electron microscopy) image of intervertebral disk cells seeded in a 3D collagen gel (gel surface)
Picture supplied by Karin Wuertz, University of Zurich, Switzerland