

Talk to us @ SLAS2013

Tecan innovation is everywhere, from state-of-the-art pharmaceutical development facilities to the smallest academic laboratories. As we see it, our job is to simplify yours, so visit Tecan on **booth #701** at SLAS2013 to see how our comprehensive range of laboratory automation, detection and OEM liquid handling solutions could benefit your laboratory.

TouchTools™ 3.0 – fully customizable touchscreen control

The latest generation of Tecan's TouchTools touchscreen user interface makes programming and operating Freedom EVO® workstations easier than ever before. TouchTools 3.0 introduces a fully customizable user interface, offering greater flexibility for instrument programmers and a structured, user-friendly environment for operators.

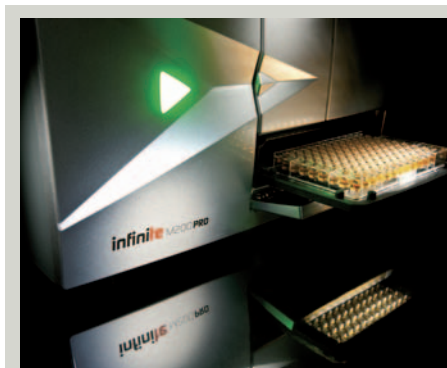


TouchTools 3.0 gives programmers the tools to create custom Maintenance and Run-Time workflows. New features make it easy to build tailored input screens, give set-up tips, and display step-by-step graphical instructions. These tools help to ensure smooth day-to-day operation with fewer operator errors. Creating assay-specific interfaces allows every aspect of a run to be described in terms familiar to laboratory scientists, including prompts, status updates, and user input screens. Embedding laboratory SOPs into the protocols will help improve compliance, enhance process security and provide traceability.

TouchTools 3.0 has a new look and feel that echoes the latest consumer touchscreen devices. This intuitive interface makes experimental set-up quick and easy, and simplified workflows reduce the time required to train new operators.

Tecan Academy – on-demand, web-based training

Going beyond traditional training courses, Tecan will soon be introducing an online training platform to provide instant access to user training and certification for various Tecan products. The Tecan Academy will offer users an easy way to learn new skills at a time and pace that suits their individual needs. Due to be launched in the US in Spring 2013, this straightforward web-based tool will enable participants to get up to speed faster, and ensures more consistent, error-free operation of Tecan hardware and software.



Infinite® 200 PRO – enhanced cell-based assays

The patent pending GCM™ (Gas Control Module) for the Infinite 200 PRO is a breakthrough for cell-based experiments, allowing simultaneous control of atmospheric O₂ and CO₂ levels within the reader's measurement chamber for the first time. Combined with the instrument's continuous shaking and heating functions, this makes the Infinite 200 PRO the first hybrid incubator/reader on the market. Recent studies have demonstrated the system's 3D assay capabilities using InSphero's organotypic microtissue model, and artefacts resulting from evaporation can be minimized by using Thermo Scientific's innovative Nunc™ Edge 96-well plate, offering a truly comprehensive solution for cell-based studies.

To find out more about these exciting new releases, visit Tecan on **booth #701** at SLAS2013