

Tecan presents new Infinite[®] 200 PRO at Forum LABO & BIOTECH 2010



Tecan chose the 9th biennial Forum LABO & BIOTECH, held at the Paris Expo Exhibition Center in Porte de Versailles, for the presentation of the new Infinite 200 PRO. The highly successful exhibition for analysis, control, research and analytical process professionals took place over four days from 1st – 4th June, and was attended by 8,564 visitors, many of them international. Tecan was one of 297 exhibitors taking the opportunity to showcase new and existing products to visitors, ranging from scientists responsible for initiating projects who were seeking information, to those who simply had a keen interest in technology.

Visitors to the Tecan stand were able to see the new Infinite 200 PRO, a system compatible with the patent-pending NanoQuant Plate for nucleic acid and protein applications. Its enhanced fluorescence bottom reading with Optimal Read (OR) function increases sensitivity for inhomogeneous adherent cells and the automated z-focus offers the highest sensitivity, even with auto-fluorescence media. Also on display was the Infinite M1000, offering flexible wavelength selection and a wide range of adjustable bandwidth

settings, high sensitivity premium Quad4 Monochromators™ technology and 3D scanning.

Liquid handling was represented by the Freedom EVO 200, with its capacity to integrate other devices on, around and even underneath the worktable. Its MultiChannel Arm™ 384 enables on-the-fly exchange between different pipetting formats for greater flexibility, and up to three arms can work independently to generate faster assay results through the ultimate parallel processing. Fast and flexible pipetting, combining eight tip pipetting across a microplate and access to individual samples with a single versatile tip was demonstrated by a Freedom EVO 75 with an 8 Plus 1 Access™ arm.

Delegates were extremely complementary about the quality of the show, the range of products represented and the contacts made there, while exhibitors appreciate the quality of visitors to this event.

