

Unique flexibility from the Freedom EVO[®] 75

Tecan has launched its latest development for the constantly evolving, compact Freedom EVO[®] 75 robotic workstation. The 8 Plus 1 Access[™] tool is a clever and cost-effective eight-channel arm for fast and flexible pipetting. This unique new arm can pipette across a microplate with eight tips simultaneously, and access an individual sample with a single versatile tip, all on the same tool.

The new option combines the best of Tecan's liquid handling experience, giving customers:

- the security of liquid level detection, down to 50 µl of conductive liquid (only for the first tip);
- the flexibility of a large volume range: 2-1000 µl;
- the efficiency of multi-pipetting with 1000 µl tips;
- the reliability and precision of the Freedom EVO series.

The 8 Plus 1 Access arm makes routine pipetting tasks, including reagent distribution, serial dilutions and plate reformatting, fast and easy to set up, and

is ideal for applications such as ELISA, PCR set-up and nucleic acid extraction. For example, fitting a Freedom EVO 75 with this new option and a robotic manipulator (RoMa) arm, offers a very economical solution for semi or full automation of ELISAs, reducing the time-consuming reagent addition steps by more than 80 % compared to using a single tip instrument. The simple and efficient Freedom EVO 75 can be equipped with on-board incubation, plate storage, plate washing and plate reading all controlled seamlessly through our easy-to-use Freedom EVOware[®] software.

The 8 Plus 1 Access further enhances the outstanding liquid handling performance, reliability and flexibility of the Freedom EVO 75, which is the perfect solution for laboratories with limited space and is ideal for those labs taking the first steps towards automation.

The 8 Plus 1 Access arm is scientific instrumentation. Not for use in human clinical or diagnostic procedures.



Capabilities of the 8 Plus 1 Access arm; eight-tip pipetting and one tip independent action

The new Infinite[®] 200 NanoQuant is launched, tailored for absorbance applications

The Infinite[®] 200 NanoQuant, Tecan's highly sensitive absorbance instrument, is being launched at LabAutomation 2008, booth 305, 27-30 January in Palm Springs, USA. The Infinite 200 NanoQuant is the first microplate reader to be developed specifically for absorbance applications with low volumes.

The new instrument represents a simplified version of the standard Infinite 200 and is available with Tecan's new NanoQuant Plate[™], a sensitive, quartz-based tool that can measure up to 16 samples simultaneously from

just 2 µl sample volumes, and detect DNA concentrations as low as 1 ng/µl. The NanoQuant Plate is compatible with multichannel pipettes and gives high measurement reproducibility; it is also available for standard Infinite 200 instruments.

The Infinite 200 NanoQuant reader is available with quad4 monochromators[™] (Infinite M200 NanoQuant) that provide high wavelength accuracy for endpoint measurements over the entire spectral range, or with new 260 and 280 nm filters (Infinite F200 NanoQuant) with a UV-



stable coating that provide exact central wavelengths and narrower bandwidths.

The instrument can perform absorbance measurements with standard microplates (6- to 384-wells) as well as half-well 96-well plates, and is easy to run with the latest i-Control[™] software that features tools for rapid DNA/RNA quantification. The Infinite 200 NanoQuant can be upgraded with the additional detection modes that are available with the standard Infinite 200 as required.