

## Tecan's Freedom EVOLyzer® becomes SFDA-compliant in China



Automating ELISAs with the Freedom EVOLyzer

Tecan has successfully registered its automated ELISA analyzer, the Freedom EVOLyzer®, with the State Food and Drug Administration (SFDA) in China. New Chinese National Standards for medical electrical equipment were introduced in May 2006 and the Freedom EVOLyzer is one of the first clinical devices to become SFDA-compliant, meaning that Tecan is leading the field in its rapid implementation of the new regulations. Seven Freedom EVOLyzer workstations were immediately ordered for customers in China following the registration. The instrument's high speed, flexibility

and walkaway features are proving to be particularly important for hospitals and blood bank laboratories, and the automated ELISA platform is currently unrivaled in China.

Mark Wang, Chief Representative of Tecan China, explained: "Tecan is well equipped to serve customers in China, with a permanent office in Beijing established in 2005, as well as a strong distribution network and nine highly trained application specialists to support the customer base."

The Freedom EVOLyzer has been designed to comply with the 98/79/EC directive on

*in vitro* diagnostic medical devices. The workstation already has plenty of satisfied customers in Europe, offering a wide range of automation options for all typical ELISA steps, including sample distribution, reagent pipetting, incubation, washing, reading and data generation. The new Freedom EVOLyzer software version 1.5, released in July 2006, includes improved error handling, more flexible scheduling and 20% faster pipetting capabilities.

*Not all of the above mentioned applications have been cleared for use in all countries, please contact your local sales office for specific information*

## Compact automation with the MultiChannel Arm 96 Gripper

Tecan's MultiChannel Arm module is now available with a Gripper as an optional add-on. The Gripper is a small module which fastens to either side of the head of the MultiChannel Arm, with fingers to grip objects such as plates and tube racks to move them around in the workstation. The Gripper can also carry out simple manipulations, eg. assembling and disassembling the Te-VacS™ vacuum separation module, giving even more efficiency and flexibility to the platform without the addition of another arm to the system.

The Gripper is an economic solution for a platform with a small to moderate throughput of 10 to 20 plates per day, and is ideal for users on a limited budget, as well as users who need a fully automated system but are reluctant to commit to a large set-up with three independent arms.

