

Tecan showcases the Freedom EVO[®] workstation configured for ELISAs to delegates visiting from Chinese blood centers

Delegates from blood centers across China recently enjoyed a visit to Tecan's European development facilities in order to share their experiences and give their input on performing high quality ELISAs within blood banking laboratories. Tecan demonstrated a Freedom EVO workstation specifically configured to meet the ELISA throughput demands observed in the Chinese clinical market, incorporating a range of features to streamline the blood testing workflow.

Efficient throughput is critical in hospital and blood center laboratories, requiring reliable automated solutions which are closely matched to laboratory workflows. The Freedom EVO offers a flexible configuration which can be tailored to individual throughput requirements, with a choice of either one or two liquid handling arms, up to two HydroFlex™ microplate strip washers, a Robotic Manipulator (RoMa) Arm for plate transportation and an Infinite[®] F50 microplate reader. The specific requirements of the scientist's assay can be fulfilled by the use of a variety of incubators, hotels and carriers, providing complete automation of the processing workflow.

The Freedom EVO is controlled by Tecan's intuitive Freedom EVOware[®] software, and uses the TouchTools Suite™ graphical user interface (GUI) to simplify day-to-day operation of the instrument. Combined with Tecan's new QC Kit for liquid handling performance verification testing, this flexible solution provides hospitals and blood centers in China with a straightforward and efficient way of automating testing.



Delegates from Chinese blood centers visit Tecan

The visit was arranged in response to increasing demand for automated laboratory systems in Chinese hospitals and blood centers, and reflects Tecan's close collaboration with customers. This year's event built on the success of the previous visit to Tecan's Männedorf headquarters in Switzerland in 2010. In addition to seeing Tecan automation in action, delegates were given the opportunity to visit Tecan manufacturing sites in both Männedorf and Salzburg, Austria, as well as a visit to a Zurich blood bank to foster relations and exchange blood bank practices. The visit also included workshops on various aspects of potential

development, providing customers and Tecan specialists with an opportunity to discuss current and future applications, as well as the bottlenecks in processing, helping to drive forward identification of novel solutions for this important market.