

Training brings the World together



At Tecan, we pride ourselves on an excellent reputation for training and, in the past year, our expertise in this area has helped to develop and encourage good relations with regulatory authorities and governing bodies all over the World.

In May 2005, Tecan was invited by China's State Food and Drug Administration (SFDA) to carry out quality system regulation (QSR) training for its members. A delegation comprising Mark Wang, Tecan's Chief Representative in China, Günter Weisshaar, the Vice President of Quality Assurance and Regulatory Affairs, and Christie Rice, Tecan's Global

Compliance Director and former employee of the US FDA, traveled to the SFDA Center for Drug Evaluation in Beijing to conduct the workshop. The course covered medical device regulations based on US guidelines, including complaint handling, adverse event reporting, corrective and preventive action, recalls and market withdrawals, all in accordance with the 21 CFR part 820 FDA regulations.

The 28 participants, all employees of the SFDA, are responsible for building up Chinese regulations to match those in the US and Europe, beginning with a Chinese Compulsory Certificate for registration of new products. The information they took from the course will help to establish national guidelines for investigating incidents when something goes wrong, whether that's

with medical devices or the companies supplying them. The workshop was very well received and was recorded and published on the SFDA homepage. Tecan also received official thanks from the SFDA and has been invited back for a second course to take place in January 2006. This workshop, for a smaller group of SFDA employees, will concentrate more closely on how to investigate complaints, the procedures involved and how to statistically evaluate the whole reporting structure.

We are delighted to have been invited back to China and will continue to use our expertise in training, whether that's in compliance and regulatory issues or in product and software workshops, to build and strengthen relations with customers, business partners and government bodies alike.

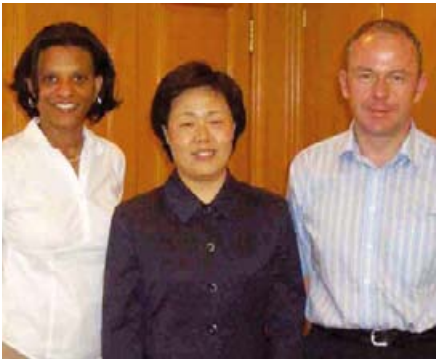


The 28 SFDA participants at the Tecan workshop
Tecan Journal 1/2006

Demonstrating the new Infinite™ and Freedom EVO® in China

CHINA

Tecan was proud to introduce Infinite, the new multimodal microplate reader, and the Freedom EVO liquid handling platform to distributors at a recent training meeting in China. The five-day event was organized by the Tecan Group Beijing Representative Office and took place in the China World Hotel (Dec 5 2005) and the Jianguo Hotel Beijing (Dec 6 to 9 2005).



Christie Rice (Director Global Compliance, Tecan US), Mrs Yi Mei (Director of Blood Management Division, Ministry of Health) and Günter Weisshaar (Executive Vice President Global Quality & Regulatory Affairs, Tecan Group Ltd.)



Mr He Ju-Cheng (General Manager of Vastec, the Tecan China distributor in the Hospital Segment), Dr Zhao Rong-Ping (Head of Immunology Division of the Clinical Diagnostic Laboratory at Shanghai Renji Hospital) and Mark Wang (Tecan Beijing Representative Office)



Mark Wang and Christie Rice

Participants came from Tecan's distributors for China, including Eastwin, Union Net Credit and Chinawealth, and Tecan's distributors for Singapore, including In vitro Technologies, Siber Hegner Sdn Bhd, SciMed Asia Pte Ltd., Everlight Biotech Co Ltd., DKSH and Pozitronics. Tecan was represented by delegates from Switzerland, Austria, Singapore, the US and India. Invitrogen was also invited to participate in the training meeting and to explore the possibility of further co-operation with Tecan.

Half of the week's meeting was devoted to the Infinite, which was introduced by Dieter Popp and Gerlinde Zerza-Schnitzhofer, and included comprehensive, hands-on training for all technical aspects of the plate reader, its software and its applications, as well as sales advice.

The product managers Jason Meredith and Christoph Jung then took the floor with the new Freedom EVO, and introduced the audience to the benefits of the new Freedom EVOware® features as well as a variety of biopharma applications appropriate to the platform.



Hands-on training demonstrations for Tecan's distributors