

Tecan enjoys success at China Forum on Transfusion Medicine 2010

Tecan recently attended the 5th National Congress of the Chinese Society of Blood Transfusion (CSBT) and China Forum on Transfusion Medicine 2010 at the Pride International Conference Centre of Chengdu in Sichuan Province, China. This international event attracted more than 1,000 experts from around the world to discuss the current challenges, trends and innovations in blood transfusion medicine. The symposium provided an excellent opportunity for both Chinese and international delegates to exchange ideas and share their experiences, and was accompanied by an exhibition showcasing the latest medical devices and technologies designed to support this sector. Tecan China worked closely with its local distributors to demonstrate the many advantages of automated liquid handling processes in this area, enjoying a steady flow of visitors to watch its state-of-the-art solutions, such as the powerful Freedom EVOLyzer® ELISA processing workstation, in action.

Alongside the main symposium and exhibition, Tecan hosted a satellite symposium on quality assurance and automated process security in the clinical environment. Günter Weisshaar, Head of Quality Assurance and Regulatory Affairs at Tecan, gave a presentation on the Company's medical device product life cycle, detailing



Tecan representatives and guests at CSBT (left to right): Yingguang Wu, Yuan Zhang, Jinming Li, Günter Weisshaar, Kirsti Tavast, Cary Shang

how this improves patient safety while better meeting the processing and regulatory needs of its customers. Tecan also invited Mr Jinming Li, Director of the Immunization Division and the National Center for Clinical Laboratory, Chinese Ministry of Health, to speak on the importance of automation during ELISA processing and genetic testing in blood banks, providing the audience with a user's perspective on the many advantages of laboratory automation.

The satellite symposium builds on Tecan's success in China as a leading supplier of sampling and ELISA processing instruments to the CSBT, and follows the recent visit to Tecan's headquarters in Männedorf by the directors and presidents of blood centers and blood banks across China to share experiences and exchange ideas.



Mr Jinming Li speaking at CSBT



Libby Zhu on the Tecan booth at CSBT