

Tecan enjoys success at 2012 SLAS Asia Conference and Exhibition

Tecan attended the 2012 SLAS Asia Conference and Exhibition at the Grand Hyatt Shanghai, China, from 19th to 21st of June 2012. This international event hosted approximately 350 delegates over three days, bringing together research scientists and business leaders from academia and industry to discuss this year's topic: *Advances in Drug and Life Sciences R&D Through Laboratory Technology*. The Conference's educational program featured renowned local and international speakers, with presentations covering a wide range of topics related to drug discovery, from high throughput screening to the use of stem cell technologies.

Tecan China demonstrated the many advantages of automated liquid handling processes in drug discovery and life science applications in the exhibition hall, enjoying a steady flow of delegates to its booth. A number of state-of-the-art automated solutions were in action on the booth, including a powerful Freedom EVO[®] drug discovery workstation equipped with a MultiChannel Arm[™] 384 (MCA 384), the top class Quad4 Monochromators[™]-based Infinite[®] M1000 PRO microplate reader and the innovative HP D300 Digital Dispenser. Tecan China's strong multidisciplinary team

of sales, service and applications specialists was joined by both Andrea Beccari – Global Head of Sales, Service and Commercial Operations – and Head of China and South East Asia Raymond Chan, and the event provided many opportunities to speak to key opinion leaders in the field in a relaxed environment.

Alongside the main conference and exhibition, Tecan hosted a tutorial on the launch of two new products. The new Infinite M1000 PRO multimode microplate reader with AlphaScreen[®] and AlphaLISA[®] technology was introduced by Marketing Application Specialist at Tecan Austria Dr Christian Oberdanner. Christian, who arrived at SLAS fresh from a highly successful series of roadshows to promote this powerful new system, explained the benefits of the system's highly acclaimed absorbance, fluorescence and luminescence scanning capabilities for academic, biotechnology and pharmaceutical research. The HP D300 Digital Dispenser was also launched at the event by Dr Yingguang Wu, Head of Marketing for the Asia Pacific region. This revolutionary sample preparation tool has the potential to streamline the drug discovery process, and received widespread interest from delegates.



Tecan staff and customers were able to discuss the many drug discovery applications of the Freedom EVO workstation



Dr Christian Oberdanner introduced the new Infinite M1000 PRO reader

The 2012 SLAS Asia was hailed a great success for Tecan, once again highlighting the Company's innovation-driven solutions and customer-focused approach to laboratory automation for drug discovery and life science laboratories.

PRODUCT NEWS

Increased flexibility and choice for Cavro[®] Centris Pump

The state-of-the-art Cavro Centris Pump is now available with a range of glass syringes and plastic valves to complement the existing portfolio of long-life ceramic components. These new options provide customers with a greater choice of materials to suit their instrument design, bringing the advanced features of the Cavro Centris Pump to a wider range of applications.

The Cavro Centris Pump is a compact, UL-recognized syringe pump module for use in OEM laboratory instruments. Offering exceptional liquid handling characteristics – with flow rates from 5 nl/s up to 5 ml/s – it sets new industry standards for accuracy,

reliability and precision. A hallmark of the Cavro Centris Pump is its elegant and robust drive mechanism, offering a broad dynamic range from a single syringe size, and the advantages of this innovative design are now available to customers wishing to use cost-effective glass syringe and plastic valve options. Already widely used throughout the Tecan Cavro range, these high quality glass and plastic components retain the premium performance benefits of the Cavro Centris Pump, while reducing the installed cost for applications which do not require the exceptional durability offered by a ceramic fluid path. Now, more than ever, you can truly tailor your Cavro Centris Pump to your instrument and application requirements.

To find out more on Tecan's Cavro Centris Pump, visit www.tecan.com/components



The Cavro Centris Pump offers a choice of glass or ceramic syringes