

AACC Clinical Lab Expo 2010

The 2010 American Association of Clinical Chemistry (AACC) Clinical Lab Expo, held at the Anaheim Convention Center in Anaheim, CA, from July 27th-29th, continues to be the world's largest show dedicated to clinical laboratory sciences. More than 21,000 people attended this year's premiere event, with an all-time record of 1,996 booths and more than 650 exhibitors.

As a leading global provider of laboratory instruments for clinical diagnostics, Tecan showcased its diversity and flexibility as your partner of choice for world class OEM and end-user solutions for laboratory automation. Highlighting this flexibility was the generic turnkey ELISA system, the Freedom EVOlyzer®, available off-the-shelf with the Tecan logo or customized with your company colors and insignia.

The Freedom EVO® liquid handling platform, combined with Luminex® Corporation's Luminex 200™ and FlexMAP 3D™ readers, provides the next dimension in multiplexing, automating xMAP® bead-based assays. Multiplex your throughput with the Freedom EVO for maximum efficiency, and automate the entire assay workflow including sample distribution, bead addition, mixing, incubation, bead washing and analysis.



Tecan Freedom EVOlyzer

New horizons in microplate detection and separation were demonstrated by the Infinite® F50 absorbance reader, a new state-of-the-art ELISA microplate reader with the latest LED technology. Also on display was the Infinite 200 PRO multimode reader, available with Tecan's NanoQuant Plate™, for absorbance applications with low sample volumes, and the HydroFlex™ 3-in-1 microplate washer, for ELISA assays, magnetic beads and cells.

Tecan's front-end automation solutions were represented by the FE500pro™. With its small footprint, the FE500pro is an effective pre-analytical automation solution which offers functions such as pre-sorting, centrifugation, volume check, clot detection, decapping, secondary tube labeling, aliquoting and destination sorting into analyzer racks, all in one system.

With Tecan OEM components, innovation was seen everywhere at the 2010 Clinical Lab Expo. A lucky attendee from Kollman Inc. correctly spotted the number of exhibitors offering products that contained Tecan components, and won an iPod®. As global leaders in pump and robotic innovation for many decades, our cutting-edge designs and



product reliability are just two reasons why brands such as LipoScience (Vantera™) and Luminex (MAGPIX®), both present at AACC 2010, trust Tecan.

Whether you're a professional in need of laboratory automation, or a partner looking for OEM solutions, Tecan provides high quality, reliable world-class solutions. We hope you'll join us and be a part of the 2011 AACC Clinical Lab Expo in Atlanta!

Luminex, xMAP and MAGPIX are registered trademarks, and Luminex 200 and FlexMAP 3D are trademarks, of Luminex Corporation.

Vantera is a trademark of LipoScience.

iPod is a trademark of Apple Inc., registered in the US and other countries.



Tecan exhibition at AACC 2010