

LabAutomation 2011

This year's ALA LabAutomation Conference celebrated 15 years of bringing together science, technology and industry to showcase the latest developments in laboratory automation. Held in Palm Springs Convention Center, California, this annual event provides a unique opportunity to meet like-minded colleagues from around the world, while exploring exciting new developments in laboratory automation. With attendance up on last year, Tecan enjoyed another successful exhibition, and customers old and new visited the Tecan booth to discover the Company's latest offerings for the life science sector.

A selection of automated workflow solutions based on the Freedom EVO® platform were in action, including automated library preparation for next generation sequencing, as well as the latest detection products, such

as the HydroSpeed™ plate washer and the patent pending Gas Control Module (GCM™) for the Infinite® 200 PRO reader. Visitors were also able to see the Attana range of quartz crystal microbalance systems on the Tecan booth for the first time. Catering to the biopharmaceutical sector, Attana's molecular characterization instruments perfectly complement Tecan's liquid handling portfolio, and users of these systems in North America will now be able to benefit from Tecan's extensive service and support network (see Corporate News on page 5).

Continuing the partnership theme, Tecan was joined on the booth by Sword Diagnostics, who introduced delegates to the innovative Sword™ Detection System, a high performance immunoassay reporter technology which takes advantage of the flexibility of Tecan's Infinite M1000 and Infinite M200 PRO series readers. Tecan instruments could also be seen on a number of other booths throughout the exhibition hall, including Artel's, where the Company's Multichannel Verification System (MVS) technology for liquid handling performance validation was being demonstrated on a Freedom EVO platform.

Alongside the main exhibition, a series of industry-sponsored technical workshops provided delegates with greater insight into the day-to-day application of these latest advances in laboratory automation. Tecan continued its support for this initiative by hosting a session on interfacing with Freedom EVOware®, as well as providing support for an MVS workshop by Artel, and we are already planning for next year's show!

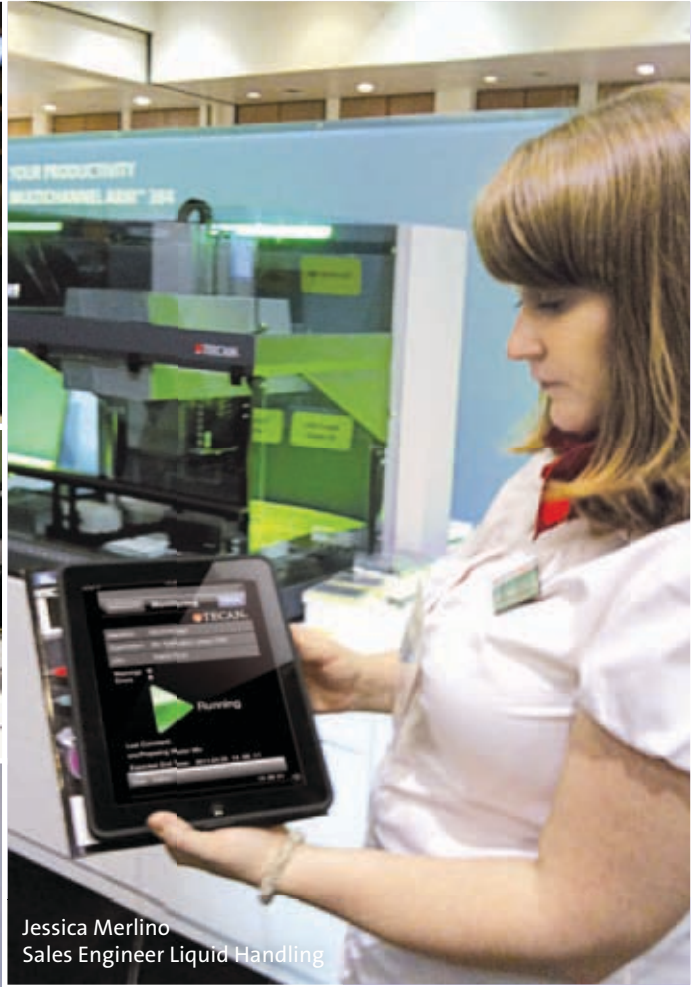




Caroline Huber
Senior Market Manager Genomics & Forensics



Read more about Tecan's partnership
with Attana on page 5



Jessica Merlino
Sales Engineer Liquid Handling

