

Tecan hits the road

Over the last few months, Tecan application specialists around the world have been organizing roadshows and seminars to discuss the benefits of laboratory automation with customers from the biotechnology, pharmaceutical and academic sectors.

In the US, Tecan has jointly hosted a series of bioprocessing seminars. Working with applications partners Atoll and FortéBio, these meetings focused on use of Tecan's Freedom EVO® liquid handling workstations in the bioprocess sector. Tecan representatives Kevin Moore and Daniel Leach presented a summary of Tecan's automation options for the bioprocessing laboratory, covering topics such as cell transfection, high throughput process development and sample preparation for HPLC, MALDI-TOF and ELISA. They were joined by Atoll's Tim Schroeder, who illustrated how MediaScout® RoboColumns can be used with the Freedom EVO platform for both process development screening and process analytical technology (PAT). Islay Campbell and Will Riches from FortéBio continued this theme, explaining the principles of the FortéBio Octet® 384 platform's BioLayer Interferometry technology, and its integration with Tecan systems.

For the Scandinavian market, Tecan's Didrik Rosen and Roland Durner hosted seminars on cell culturing in Copenhagen and Stockholm. Invited speakers from leading companies in the field shared their experiences of combining experimental theory with practical applications in the laboratory, introducing attendees to the latest developments in cell culturing technology. There were also workshops to demonstrate the practical advantages of automation for a range of applications, illustrating how easy it can be to implement laboratory automation, even without previous experience, using Tecan's innovative TouchTools Suite™ touchscreen control system.

Finally, representatives from Tecan's UK offices have been touring England and Scotland in a series of one day exhibitions. This roadshow introduced the latest laboratory automation technologies and multifunctional microplate reader designs to a broad scientific audience, visiting universities, research institutes and science parks. Visitors were able to discuss the huge wealth of applications which can now be automated, and the potential benefits of Tecan's technology, including increased throughput, improved reproducibility and greater confidence in results.

To find out which events Tecan will be attending in the coming months, see the back cover or visit www.tecan.com/events

MediaScout is a registered trademark of Atoll GmbH, Weingarten, Germany.

Octet is a registered trademark of FortéBio Inc.

