

Tecan UK customers gather at historic Hinxtton Hall

The Tecan UK User Meeting 2013 provided the perfect opportunity for members of the automation community to network with other users, discussing a wide array of applications.

Towards the end of 2013, Tecan UK hosted a two-day user meeting at the Wellcome Trust Conference Centre, home to the world-renowned Sanger Institute, in Hinxtton. Delegates at the state-of-the-art venue, located in over 100 acres of parkland on the banks of the River Cam, were able to enjoy the facilities of the purpose-built conference center and the historic Hinxtton Hall. Held from the 5th to the 6th of November, the event provided the opportunity to hear speakers from many different fields – including the pharmaceutical, biotechnology and academic sectors – discussing the

wide variety of applications for which Tecan platforms are used, complemented by poster presentations and a display of some of Tecan's automated solutions.

Presentations covered the adoption of automation in diverse areas such as lab-scale protein production, high throughput ELISA, biobanking and protein-DNA interaction screening, as well as the challenge of incorporating sample collections obtained during company acquisitions into one standard format. A highly informative presentation on the theory behind optimizing liquid classes – supported by a presentation on the practicalities of implementation in the laboratory – was well received, and there were tips on maintaining Freedom[®] EVO systems and using Freedom EVOware[®]. Delegates also heard about the role of the Tecan Integration Group in providing customized solutions that are a little outside of the ordinary, as well as new Tecan applications. The event concluded

with a lively round table discussion, where customers and Tecan specialists debated the use of automation, its challenges and the solutions.

The success of the event was summarized by Steve Hawthorn, Commercial Manager, Tecan UK Ltd, who said: "User meetings are about supporting our customers, bringing together the different branches of the life science market and creating an automation community where our customers can network and learn from each other's experiences. The varied program attracted scientists from a broad range of backgrounds, and there was ample time for networking during the day and at the evening dinner. Feedback was very positive; delegates will be recommending the event to colleagues and attending again themselves, and we have already been approached by a number of customers interested in presenting at future user meetings."

"Feedback was very positive; delegates will be recommending the event to colleagues and attending again themselves."

