

# Systems biology symposium, a great success

Our systems biology symposium, held in mid-September on the shores of Lake Zürich, presented a rare opportunity for leading scientists working at the forefront of research to meet together in an informal atmosphere. The event was attended by researchers representing many of the world's major research institutes, and several prominent scientists gave presentations on their current and future initiatives in this fast-moving field.

The event was arranged in association with Alan Sawyer, Head of the EMBL Monoclonal Antibodies Core Facility, who chaired the first day. He began proceedings with a presentation about

the new monoclonal antibody production facility currently under construction at Tecan's headquarters in Männedorf and on site at Monash Antibody Technologies Facility (MATF) at Monash University, Australia, where he is soon to take up a position as Director.

Alan said: "The timing of this event was perfect. Some colleagues and I had discussed how we could bring together top class scientists working in systems biology and, almost simultaneously, Tecan approached me expressing an interest in just such a forum. The result has been an intensely stimulating environment, with plenty of opportunity for in-depth, constructive discussions on a more

intimate level than is normally possible at larger scientific meetings."

Our Senior Vice President, Carl Severinghaus, said: "The entire event has been a very rewarding experience and it was a privilege to bring so many luminaries together and hear first hand about their latest projects. We have also been able to learn more about the role of automation in systems biology and have a better understanding of how Tecan can provide the ideal automation solutions for these applications."

A similar event is already being planned for 2008.



*Participants at the Tecan Symposium 2007*