

Tecan symposium 2008

'Biologics: from benchtop to production'

Tecan is keenly looking forward to hosting the second Tecan symposium on September 16-19, 2008 in Oerlikon, Zürich, Switzerland, following last year's hugely successful inaugural event. The annual scientific forum offers an intimate opportunity for leading scientists and a varied range of Tecan users to meet and discuss emerging scientific issues in connection with Tecan's fields of expertise.

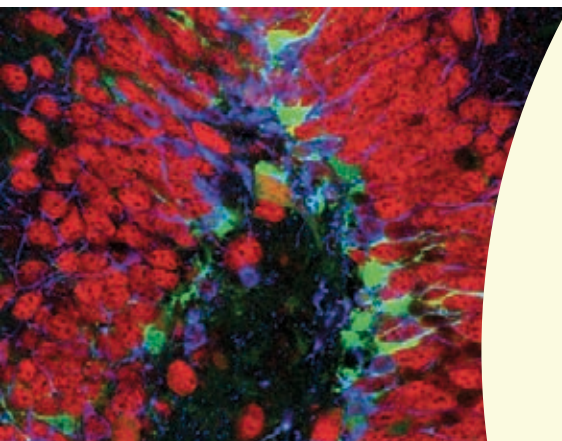
The 2008 symposium, entitled: 'Biologics: from benchtop to production', will cover the rapidly growing areas of biotherapeutics, stem cell research and regenerative medicine. The formal program will be held over four sessions:

- advances and opportunities in stem cells
- from cells to therapy
- high throughput applications
- optimization and production of therapeutic proteins and antibodies.

Scientists at the forefront of their respective fields will present on various cutting-edge initiatives in research and application of biologics, followed by open, in-depth discussions with the invited audience.

The first Tecan symposium was held in September 2007 in Erlenbach, Zürich, Switzerland, focusing on the subject of systems biology, and was chaired and jointly organized by Alan Sawyer, director of the Monash Antibody Technologies Facility in Australia. Alan will return in 2008 to present his valued input on high throughput monoclonal antibody production. With many luminaries in research and industry already confirmed to attend, this year's meeting promises to be another successful event, and is eagerly anticipated by all at Tecan.

www.tecan.com/symposium2008



The different cell populations in the hippocampus of the adult mouse during neurogenesis. Image taken with a confocal microscope at 40x magnification. Image supplied courtesy of Professor Kempermann, CRTD.

Zürich **2008**
second Tecan
Symposium