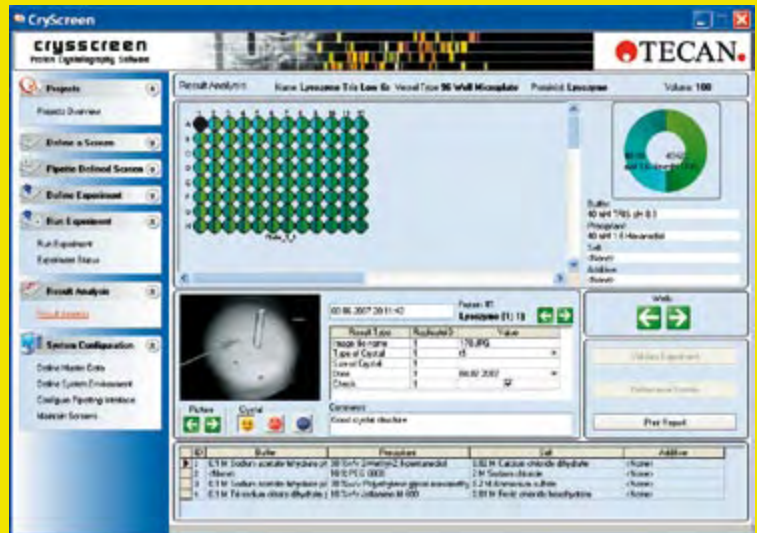


Enhance your crystallography with the new CrysScreen™ 2.0 software



An upgraded version of the software offers a number of user-friendly improvements in the management of protein crystallography trials.

Tecan has recently upgraded its CrysScreen protein crystallography software. The new version 2.0 runs under Tecan's Freedom EVOware® standard software, ensuring simple and efficient crystallography process control using the Freedom EVO® liquid handling workstations. It contains a number of user-friendly improvements, including a data export function. The CrysScreen package provides easy management of protein crystallography trials, including buffer creation, set-up of experimental conditions, running of experiments, data management and analysis of results. Watch out for further information on CrysScreen 2.0 very soon!



Work starts on the largest monoclonal antibody production facility in the Southern hemisphere

Tecan and the Monash Antibody Technologies Facility (MATF) at Monash University, Australia, have announced a partnership to create one of the largest, most sophisticated monoclonal antibody production systems in the world. Plans are already underway to develop the system, which will be based around multiple, integrated liquid handling workstations and will fully automate all the stages of the production process, as well as incorporating state-of-the-art, third party devices.

The MATF is one of the only high throughput production facilities in the world offering custom-made, high-affinity monoclonal antibodies. The new system will increase its production capacity from original estimates of 500 panels of novel antibodies a year to potentially 5,000, helping to alleviate the bottleneck of antibody supply experienced by many researchers working on protein function and physiological processes. The project is partially funded by the Victorian State

Government, the Australian Stem Cell Centre and Monash University.

Tecan's team of experienced application specialists and engineers will build the new system at the company's headquarters in Männedorf, Switzerland, ready for installation at the MATF in April 2008.



MATF workers (l to r) Director Alan Sawyer and Deputy Director Michael Spiegel