



# The 6th International BioForum & BioExpo Japan



*Dr Titus Kretzschmar (amaxes) and Eric Willmann (Tecan) presented a seminar on automated cell culture and transfection screening*

**Proclaimed as Asia's largest bio-event,** the 6th International BioForum & BioExpo Japan was held at the Tokyo International Exhibition Center in Tokyo Bay, also known as Tokyo Big Sight, on June 20-22, 2007. During the three-day exhibition there were almost 19,000 visitors, consisting mainly of researchers from universities, research centers and medicine/food manufacturers.

The International BioExpo attracted around 600 national and international companies, ranging from instrument and chemical manufacturers to suppliers of bioinformatics solutions and bioventures, supporting various life science research areas like pharmaceuticals, gene therapy, genomics and proteomics.

With 'automate your cell-based assays with Tecan' as the theme, Tecan Japan demonstrated cell-based automation solutions based on the Freedom EVO® workstation, together with other product line-ups including the latest microplate readers and washers like the Infinite™ F500 and the HydroFlex™. In the joint booth, REMP showed the Sample Safe™,

the newly released, fully automated -80 °C sample storage and retrieval system. Tecan's staff were so inundated with visitors to the booth, that some needed to book appointments in advance for the product specialists to be available for more detailed discussion on specific products and solutions for the customers' application needs. A Freedom EVO 200, fully equipped with a liquid handling arm, MultiChannel Arm™ and robotic handling arm, operated continuously during the exhibition, and the Te-Flipper™ undoubtedly attracted a lot of attention from passing visitors.

The BioForum was held concurrently with the BioExpo, with a total of 250 presentations, including keynote talks from eminent life scientists. There were also research presentations by universities and public institutes in the BioAcademic Forum, technology presentations in the BioVenture Forum, and presentations about products/technologies that support bioresearch by various exhibiting companies. Dr Titus Kretzschmar, Vice President of Research and Development at amaxes AG and Eric

Willmann, Marketing Manager, Cell and Protein Science at Tecan, made a joint presentation entitled 'Automated cell culture and transfection screening' in the Exhibitors' Product/Technology Seminar. This seminar introduced Tecan's Freedom EVO integrated with amaxes's Nucleofector® 96-well Shuttle® transfection system, and detailed automation solutions for high throughput screening and RNAi screening, fully automated cell culture, preparing cell lines, cell function analysis and cell-based assays. Many pharmaceutical companies participated and, even after the seminar, they were enthusiastically asking questions at the exhibition booth.

Nucleofector and 96-well Shuttle are covered by patent and/or patent-pending rights owned by amaxes

