

The International Union of Biochemistry and Molecular Biology (IUBMB) 2006

Tecan shows new products in Japan

The 20th IUBMB International Congress of Biochemistry and Molecular Biology, covering all topics within the fields of molecular biology and biochemistry, was held from June 18-23 at the Kyoto International Conference Hall, Takaragaike Prince Hotel, in Japan. Run once every three years in various locations around the world, it has been 39 years since the congress last visited Japan. At the opening ceremony, the 8,000 participants from 71 countries were graced by the presence of His Imperial Highness the Crown Prince.

As a joint event with the symposia of the Japanese Biochemical Society and the Molecular Biology Society of Japan, normally held in autumn, the program with the theme "Life - Molecular Integration & Biological Diversity" was aimed at giving young researchers, who carry the future of Japan's life science, the opportunity to directly experience some of the world's top class research. Prominent researchers presented 11 plenary lectures, symposia across 89 themes, and 6,000 posters.

In the Scientific Exhibition, with 271 participating companies, Tecan Japan displayed the Freedom EVO® with the



new 96-channel pipetting head (MultiChannel Arm, MCA), Cellerity™'s Flask Flipper™ module, and the Infinite™ 200 microplate reader. Nicolas Sandoz, Product Manager of the MCA, and Roland Durner, Tecan's Global Product Manager, attended the event to promote the Cellerity. After visiting Tecan users throughout Japan during the week prior to the exhibition, Nicolas Sandoz and Roland Durner offered further expertise at the Tecan exhibition booth, demonstrating products seen in Japan for the first time. REMP's Small-Size Store™ (SSS) and devices related to the REMP Tube Technology™ were also exhibited.

Many existing users of Tecan's automated systems visited our IUBMB exhibition stand to find out more about the MCA, a new option on the Freedom EVO. The Infinite microplate reader, which has been very popular since its launch last year, has maintained a high level of interest, and the exhibition offered many opportunities to obtain feedback from experienced users of the instrument. The knocking sound from Cellerity's Flask Flipper module dislodging adherent cells from the inner surface of flasks attracted much attention, prompting many people to listen to the descriptions from Tecan staff.

