

Tecan teams up with high profile business partners in a number of new developments

In recent months, even more collaborations with industry partners have come to fruition, bringing application-specific, automated laboratory systems to researchers working in a range of disciplines.



Tecan and Applied Biosystems have jointly developed the HID EVOLUTION™ System for the automated analysis of forensic DNA casework samples. This new standardized system provides fully automated sample quantitation, normalization and short tandem repeat (STR)-PCR set-up, by integrating Tecan's Freedom EVO® 150 liquid handling workstation with the Applied Biosystems 7500 Real-Time PCR System, 3130x1 Genetic Analyzer, and Quantifiler® and AmpFℓSTR® DNA testing kits. The HID EVOLUTION is expected to reduce the amount of time required to process forensic DNA samples and help laboratories to increase their sample throughputs.

Tecan and amaxa, Inc. have successfully integrated amaxa's Nucleofector® 96-well Shuttle® System with the Tecan Freedom EVO liquid handling workstation, allowing fully automated, high throughput transfection of

primary cells and difficult-to-transfect cell lines, including non-dividing neurons and T cells, with siRNA, DNA and other substrates. The development is ideal for large-scale studies that involve high throughput transfection stages, such as RNAi-based screening for drug development. The 96-well Shuttle® relies on unique electroporation technology for transfection of primary cells with high efficiency and excellent survival rates. Integrating this system with the Freedom EVO workstation entirely removes all manual steps for cell preparation and transfection, allowing 96-well microplates to be completely processed in just minutes, depending on the protocol used. You can read more about the collaboration on pages 9-11.

Tecan and VWR International, Inc. have announced a strategic alliance to unite the strengths of each company and together provide Tecan's cutting edge

microarray instruments and microplate readers to the global research market. The collaboration takes advantage of VWR's BioSciences program, which aims to provide researchers with the highest quality services and products available, whatever their application, and a comprehensive sales network – including more than 120 sales representatives worldwide and 15 BioSciences managers in Europe alone. Tecan's range of technically excellent microarray products and microplate readers offers perfect solutions for customers of the BioSciences program working in all areas of drug discovery and life science.

Tecan and Cerionx, Inc. have made an exclusive distribution agreement for the supply and marketing of the Cerionx 8-channel TipCharger™ System to the worldwide biopharmaceutical market. Addition of the TipCharger™ System as an integrated and supported component to Tecan's automated liquid handling and robotics workstations offers uniquely superior pipette tip cleaning and sterilization. The TipCharger™ System uses room temperature ('cold') atmospheric pressure plasma to clean and sterilize pipette tips by removing contaminants at the molecular level without using hazardous solvents. This revolutionary plasma cleaning technology reduces the potential for contamination at the sample and consumable levels, reduces run time and improves assay productivity. Find out more about the Cerionx TipCharger™ System on pages 20-21.

Quantifiler and AmpFℓSTR are registered trademarks of Applied Biosystems or its subsidiaries in the US and/or certain other countries.

The Nucleofector 96-well Shuttle System is covered by patent and/or patent-pending rights owned by amaxa.

TipCharger is a trademark of Cerionx, Inc.