

A new generation of Tecan OEM components

Visitors to LabAutomation will be able to see Tecan's new generation of liquid handling technology components, the new pump and the new robot.

New pump

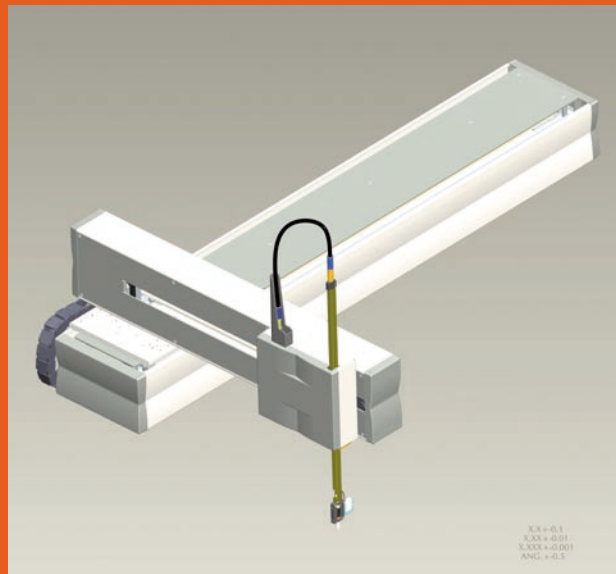
The new pump was developed to complement and extend our current range of syringe pumps, including the Tecan Cavro® XCalibur and Tecan Cavro XLP. The new pump increases standards for accuracy and precision through a new, simple but rigid direct drive pumping mechanism. The use of long-life ceramics for the fluidic path ensures unbeatable longevity, even for solvents. It also offers a broader working volume range from a single syringe size compared to other similar pumps, enabling users to do more with one pump than ever before. The new pump features increased on-board diagnostics, so the user can interrogate the pump for its critical performance criteria.

New robot

The innovative new robot provides users with the security of knowing exactly what the product is doing and where the pipetting probe is located at all times, minimizing probe positioning error. Its simple design promises the same robustness and reliability that is characteristic of Tecan's products, while its modularity offers superior flexibility, making the new robot easily configurable for many different applications. The new robot is designed for easy integration, being both easy to install and to operate, making it the ideal OEM robotics component choice. It is encased so that all working parts are hidden to sport a finished professional appearance.



New pump shown with 1.25ml syringe



New robot shown in single arm configuration