

Construct your own robot – custom design for the Freedom EVO® 75

Now you too can build your own robot. Tecan has just introduced the Freedom EVO configurator, a web-based application which enables the user to custom design a Freedom EVO 75 platform to meet a laboratory's specific requirements.

This easy-to-use website allows scientists to build a virtual Freedom EVO 75 platform by adding individual modules, such as liquid handling arms, plate handling robotics and labware carriers, to meet their exact requirements. Alternatively, the user can choose from a number of pre-defined layouts – PCR, ELISA, xMAP assay, and nucleic acid extraction – which can be customized by the addition or removal of components,

as required. A drop-down overview of the selected options enables the user to review the platform configuration and to make amendments if necessary. Once the final platform layout has been determined, the chosen configuration can be printed for future reference.

To discuss the custom built Freedom EVO 75 in further detail, or to place an order, contact the Tecan sales team via the 'Contact Sales' button on the website.

To find out more, and to design your own Freedom EVO 75, visit www.tecan.com/freedomevo75



Freedom EVO configurator allows you to construct a virtual Freedom EVO 75.

TouchTools Suite™ generates a buzz with conference attendees

TouchTools Suite has proven to be a star attraction among visitors at the numerous conferences and exhibitions Tecan has attended since its launch last year. The intuitive, touchscreen control system has made automated liquid handling more accessible to many scientists, allowing real-time control of Freedom EVO® platforms without the need to program pre-defined scripts.

Eskil Trollhagen, Tecan product manager for the TouchTools Suite, recently received a SelectScience.tv Web Award for his video guide through the system's capabilities, and explained the reasons for its success: "TouchTools Suite graphically guides the user through setting up and running a method on the Freedom EVO workstation, making it easy for inexperienced users to operate the

platform, while minimizing the risk of errors. This user-friendliness has been very well received by occasional users and automation specialists alike. For specialists it means that they do not need to guide colleagues through running even very basic methods and, for inexperienced users, it allows them to take advantage of the Freedom EVO's automation capabilities without having to invest time in automation training."

"One of the most attractive features of TouchTools is Instant Pipetting™, which allows the operator to control the platform in real time, an industry first. This has been a real draw for conference attendees, allowing them to see the Freedom EVO performing their commands immediately. TouchTools Suite is so easy to use that many visitors were actually teaching each other to use it."

For details of events where you can meet Tecan in the coming months, see the back page or go to www.tecan.com/events.

To find out more about Tecan's TouchTools Suite, visit www.tecan.com/touchtools

