

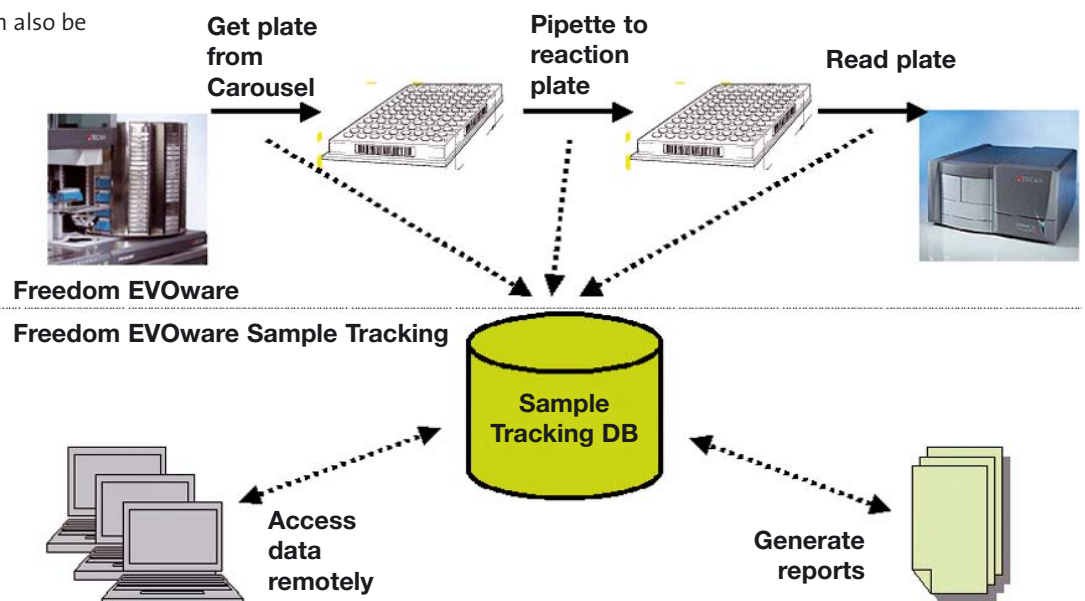
Software focus: four new packages from Tecan that unite to integrate your data

At Tecan we realize that communication between your instruments is key, so we have released a range of new software packages that enhance connectivity, for simple integration of your plate readers with Tecan's Freedom EVO® workstations and seamless combination of your data. The four new software packages are Freedom EVOware® Sample Tracking, Freedom EVOware 1.2, i-control™ and Magellan™ 6.o.

Our Freedom EVOware Sample Tracking makes it easy to electronically record the status of all your samples throughout your entire application. Sample Tracking can register scanned labware then integrate and process all data acquired for each sample, including pipetting and plate reader data. These are stored and then used to generate and export reports containing the results. Sample Tracking is compatible with an extensive range of Tecan products, simplifying the logistics of handling large numbers of samples and improving your quality controls. It also integrates data from Freedom EVOware 1.2 and Magellan 6.o for simple combination of plate readers and pipetting platforms. Sample Tracking is very much central to the interactions of all the new software (see figure), since it tracks movement of all plates (as controlled by Freedom EVOware) and all pipetting actions. Data and results from plate readings are fed to Sample Tracking for the generation of reports at the end of the process and these can also be accessed remotely.

Freedom EVOware 1.2 integrates the new i-control software (see below), for straightforward combination of the Infinite™ 200 multimode microplate reader with your Freedom EVO, and is also compatible with the new Magellan 6.o, allowing integration of other multimode plate readers with the workstation. The integration means that the Infinite reader, for example, will appear in Freedom EVOware's control bar and script editor in the same way that other supported devices appear and are called up, for easy process definition. Freedom EVOware's functionality can be extended with add-ons such as Sample Tracking, Normalization and Freedom EVOware Sample Oriented software. It supports a variety of hardware, both from Tecan and from third parties, bringing greater flexibility to your laboratory.

Freedom EVOware 1.2 has many other new, user friendly features to help your applications run more smoothly. These include a start-up wizard, which provides simple access to common daily routines, helping you to get your processes up and running quickly and easily, and the intuitive graphical user interface, which makes new assay protocols easy to create. There are also wizards to help you create advanced pipetting procedures, including relatively complex routines, and robotic moves are graphically defined, helping you to visualize processes. Sequences can be simulated and viewed, making it quicker and easier to develop new protocols as well as allowing remote monitoring. The package also comes with an expanded library of device drives and its modular architecture makes it easy to extend the software as new devices become available.





Freedom EVOware supports a wide range of Tecan modules, including the LiHa arm and dual LiHa arm, RoMa arms, Pick and Place (PnP), Te-MO™, Te-MagS™, incubators, hotels, PosID™ 2 and PosID 3. The new Freedom EVOware 1.2 additionally supports Infinite (through i-control), Sample Tracking, Sample Oriented.

Freedom EVOware 1.2 is available in two versions, Freedom EVOware Standard and Freedom EVOware Plus, which has extra benefits including a fully integrated pipetting scheduler, which can scale up your application for multiple plate runs without requiring new software. There is a dynamic scheduler for efficient use of resources and an autocalibration of schedule feature, which uses real time to create a better sample processing schedule, allowing higher throughput. The Freedom EVOware Standard allows synchronous or asynchronous pipetting with two LiHa arms, while in Freedom EVOware Plus the two LiHa arms can either be used for one process or different LiHa arms can be used for different processes.

Tecan's new i-control software installs itself into the Freedom EVOware 1.2, so that the Infinite 200 plate reader can be fully combined with other application processes running on the Freedom EVO. Typically, when there is a plate reader integrated into a Freedom EVO platform, a RoMa arm transports the sample plate from the workstation and puts it into the reader, where the plate is read and data are generated; the RoMa arm then removes the plate. Fully integrating the reader's control software into Freedom EVOware allows complete automation of the reader processes by your Freedom EVO and the i-control takes over from Freedom EVOware once the microplate is transferred into the reader. It then passes the control back when the reading is complete, so this compatibility between the software is a really beneficial new feature. The new i-control has the same look and feel as Freedom EVOware 1.2.

The new Magellan version 6.0 is compatible with all of Tecan's multimode plate readers and has powerful data reduction capabilities, including being able to perform multiplate analyses, comprehensive spectra calculations and 3D scans. It supports all major detection modes, as well as the injector system, and is designed to allow customers to comply with 21 CFR part 11 regulation, for electronic records and electronic signatures. The new program has an intuitive, wizard-guided user interface

that maximizes its ease of use, allowing seamless process control, automated data import functions in variable formats, and easy and fully automated data export to LIMS. The program even comes with example files of typical applications to get you started quickly.

For more information on our new software, please contact Jason Meredith, Global Product Manager Software: jason.meredith@tecan.com

Freedom EVOware® on Freedom EVO® Clinical, Tecan's automated platform for clinical diagnostics

The Freedom EVO Clinical with Freedom EVOware and its new module for sample oriented processing, which is predominantly used for clinical applications, is a powerful tool for controlling the Freedom EVO Clinical family of application-oriented open platforms for clinical diagnostics and blood bank environments.

The Freedom EVO Clinical/EVOware system offers:

- application-oriented, scalable precision liquid handling and robotic platform for clinical applications like EIA preparation and pooling
- extended functional spectrum by integration of new options for clinical use
- enhanced process control and safety features without compromising instrument flexibility
- scheduler function for increased throughput
- true walk-away capability

Freedom EVO Clinical is not available in the USA although sample oriented Freedom EVOware is available with pooling functions disabled.



For more information on Freedom EVO Clinical please contact Christoph Beck, Product Manager Corporate Account & Clinical Diagnostics: christoph.beck@tecan.com