

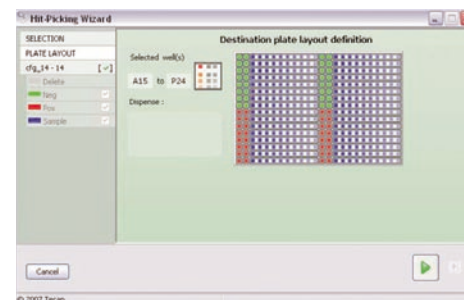
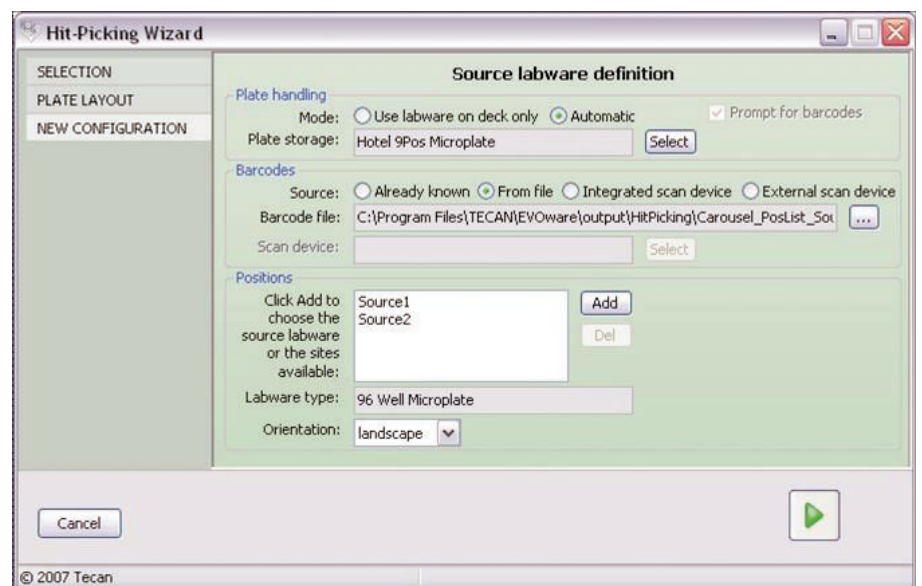
Automated cherry picking with the Freedom EVOware® Hit Picking Wizard

Tecan's Freedom EVOware® Hit Picking Wizard is available as an add-on to your Freedom EVOware software for automated plate and compound management during hit picking. The software is suitable for a variety of hit picking applications with the Freedom EVO® platform, such as production of IC₅₀ plates for hit confirmation, cherry picking after high throughput screening, or providing secondary screening plates for use in ADME assays.

Source plates and empty plates for hit compounds are loaded into storage devices or placed on the Freedom EVO deck and compounds of interest, identified in a hit list provided by the LIMS, are pipetted into the assay plates by the Freedom EVO liquid handling arm. The workstation can also prepare controls and dilutions for further assays if required, and the software provides a full report at the end of the process.

The Hit Picking Wizard integrates fully with the main graphical user interface of both the Freedom EVOware Standard and Freedom EVOware Plus programs. The software handles a variety of storage devices (eg. carousels, stackers and hotels) and labware (eg. REMP tubes and delivery plates, 96- and 384-well plates), and provides a simple graphical layout of the plate for straightforward hit picking. It controls a single 1-, 2-, 4- or 8-tip liquid handling arm, and is compatible with 200 and 1,000 µl disposable tips.

The Hit Picking Wizard also takes care of plate logistics, such as instructing a robotic manipulator (RoMa) arm to replace source plates and swap destination plates as necessary during the process. The add-on integrates with the Freedom EVOware Sample Tracking software, so that sample tracked labware continue to be identified and tracked during the hit picking processes.



Screenshots from the Freedom EVOware Hit Picking Wizard