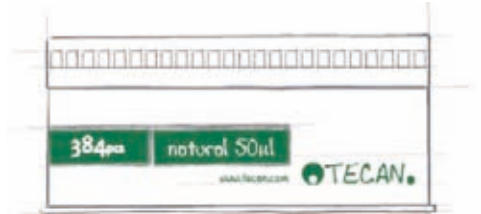


Quality disposable tips for reliable 384- and 96-channel pipetting

Tecan has introduced the new MCA384 disposable tips, complementing the recently launched versatile MultiChannel Arm™ 384 head. The system brings higher productivity to automated liquid handling applications, with reliable 8-, 12-, 16-, 24-, 96- and 384-channel pipetting in a broad volume range of 0.5 to 125 µl.

MCA384 disposable tips are available in two volume sizes – 50 µl and 125 µl. The 15 µl tip will be available in the near future. The tips have been engineered with a unique sealing system, where each tip has its own internal, built-in seal.

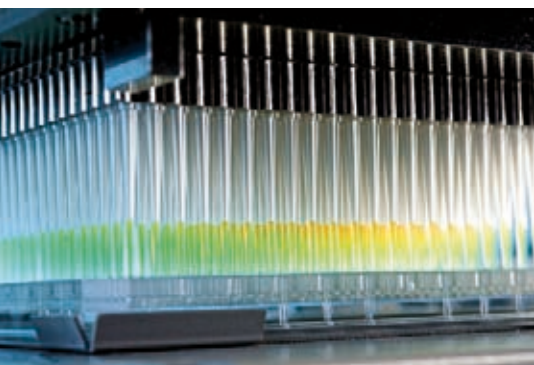
Tecan's MCA384 disposable tips are carefully designed and produced for optimal performance with the MultiChannel Arm 384 head. The tips are supplied in newly designed racks that are color-coded and clearly labeled to indicate the tip volumes for straightforward identification on the worktable. The rack labels also document lot numbers, for easy traceability.



The new rack design for easy traceability on the worktable

The flexible MultiChannel Arm 384 head is able to switch quickly between 384- and 96-channel pipetting, as well as pipetting using one row or one column of tips, even during a run, depending on the requirements of the application. The precision-built 15 µl tips allow pipetting into 1,536-well microtiter plates, while the 125 µl tips can access the bottom of a deep-well plate.

For more information about Tecan's disposable tips, visit www.tecan.com/disposablestips



Tips in action – 384 15 µl disposable tips dispensing liquid into a 1,536 microtiter plate

Freedom EVOLution™ 2.1 – sample preparation and splitting on one platform

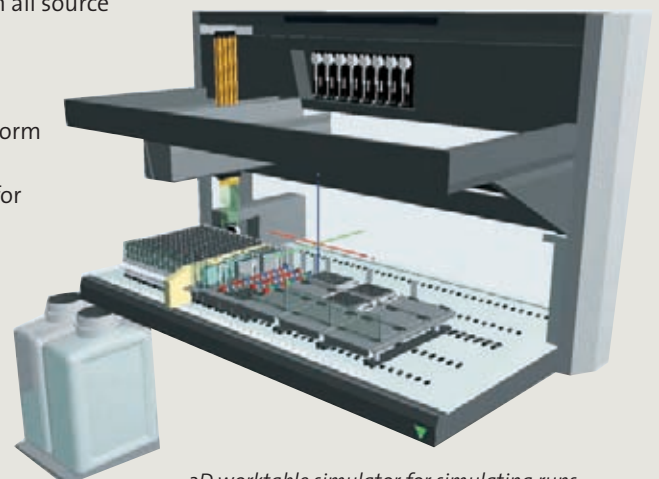
Freedom EVOLution 2.1 is the latest version of Tecan's software for Freedom EVO® Clinical platforms, offering additional application areas – sample splitting and archiving – in addition to all the existing sample preparation functions. The software allows laboratories to integrate several applications on one instrument.

The new software makes it possible to automate the splitting of source tubes into multiple destination tubes, and the archiving of source tubes into archive plates, without the need for a LIS/LIMS connection. Sample traceability is maintained by barcodes on all source and destination tubes, and the workflows are standardized, making the Freedom EVO Clinical platform ideal for all laboratories handling clinical samples for analysis or archiving.

Freedom EVOLution 2.1 is available as standard on new instruments, and as an upgrade option for existing instruments. For more information about Tecan's Freedom EVOLution, visit www.tecan.com/evolution



User actions are guided by wizard-driven GUI



3D worktable simulator for simulating runs