

Freedom EVO[®] Clinical 75 now available for nucleic acid testing



Handheld barcode scanner option

Tecan has launched the Freedom EVO Clinical 75 NAT workstation, a compact pipetting platform specially configured for extracting genomic DNA, viral DNA or RNA from clinical samples. It automates nucleic acid purification from serum or plasma, and allows flexible input from one to 24 samples; the workstation is fully compatible with the Abbott *mSample* Preparation System RNA and DNA kits. The extracted RNA or DNA showed excellent assay performance with Abbott RealTime[®] HIV, RealTime HCV and RealTime HBV viral load assays during customer evaluation. This latest addition

to the Freedom EVO Clinical family has been designed and built to meet the IVD-D (98/79/EC) requirements, as with the larger Freedom EVO Clinical 100, 150 and 200 platforms.

The instrument is equipped with a 2-tip liquid handling (LiHa) arm and the Te-MagS[™] magnetic bead separation module, with an adapted 24-position block allowing simultaneous pipetting of two samples; a barcode scanner can also be integrated with the system for sample identification. Tecan's Freedom EVOware[®] Standard software controls the instrument and allows dedicated and open-mode scripts to be run. The integrated Sample Tracking software wizard provides complete tracking and documentation of all samples throughout the process.

Not for sale in the USA and Canada.
RealTime is a trademark of Abbott Laboratories.



Optimized layout for 24 samples

Quality and quantity control for pooling of clinical samples

Tecan has recently completed its first installation of a Freedom EVO Clinical liquid handling workstation combined with the Te-PoolSafe[™] option and the Pressure Monitored Pipetting (PMP) option, controlled by Freedom EVOware Sample Oriented software. The integration of both options with the platform brings a new level of safety and quality assurance to automated blood pooling laboratories. For the first time, precise monitoring of both the dispensed volume and quality of each pipetted sample can be simultaneously performed, providing greater process security.

The Te-PoolSafe allows a fast and sensitive liquid arrival check by monitoring each pipetting step and comparing every

single dispense made with the expected volumes. The PMP tool analyzes pipetting quality by detecting faults such as clots, air aspiration, non-dispensed samples or system leakage; faults are detected by comparing recorded and real-time simulated pipetting pressure signals. The pressure simulation relies on fluidics, user-defined parameters (such as sample type and pipetting volume) and dynamic fitting of sample type and system-specific variations (eg. in sample viscosity or system liquid properties). Tecan's extensive prevalidation work of the two mainly used liquid classes - serum and water - allows quick and easy PMP installation and application set-up.

Not for sale in the USA and Canada.



Typical pooling set-up with Te-PoolSafe