

New TouchTools™ PCR Wizard aids automated PCR set-up

Tecan has launched the TouchTools PCR Wizard, an easy-to-use Freedom EVOware® add-on offering straightforward automation of PCR reaction set-up on Freedom EVO® platforms. This new tool simplifies the set-up of applications – including end-point, real-time and multiplex PCR, sequencing, genotyping and gene expression, as well as pathogen and mutagenesis detection – helping to reduce training time and costs.

The PCR Wizard's effortless touchscreen operation and step-by-step instructions guide users through the set-up process, allowing sample information to be entered manually, imported from a file or identified by barcode scanning. Users can quickly and easily optimize 96- and 384-well PCR plate layouts to meet specific workflow needs, using up to 96 master mixes and defining their own controls and dilution curves. Seamless data transfer to Bio-Rad and Life Technologies real-time cyclers avoids transcription errors,

and optional integrated sample tracking further increases process security. With just a single set-up required for multiple PCR plates per run, efficiency is considerably enhanced and, for maximum flexibility, the PCR Wizard can be used in conjunction with the Freedom EVOware Normalization Wizard.

To find out more on Tecan's TouchTools PCR Wizard, visit www.tecan.com/pcrwizard



The PCR Wizard offers easy automation of PCR reaction set-up

Nested disposable tips now available for Liquid Handling Arms



The nested format allows up to five trays of tips to be stacked on each SLAS-format carrier

The demand for ever-increasing throughput means that many laboratories are looking for solutions to increase the throughput of their automated systems. Tecan has created Nested LiHa Disposable Tips specifically to meet this need, allowing up to five trays of the new 350 µl tips to be stacked on a single SLAS-format carrier, with fully automated tip and rack handling through Freedom EVOware.

This cost-effective, space-saving solution increases labware capacity and allows optimized use of deck space, minimizing the need for manual intervention and increasing walkaway time for Freedom EVO liquid handling workstations. The nested

350 µl tip is available in both Tecan Sterile and Tecan Pure options for standard applications, with a Te-Stack version available soon in Tecan Standard purity for higher throughput applications.

To find out more on Tecan's Nested LiHa Disposable Tips, go to www.tecan.com/consumables