

Cavro[®] ADP – a new innovation for liquid handling

Tecan has launched the new Cavro Air Displacement Pipettor (ADP), a compact, pneumatic pipetting solution for easy integration and exceptional liquid handling performance. The Cavro ADP is a fully programmable, pneumatic pipetting module designed for integration into instruments requiring minimal maintenance or where space is at a premium. Using pneumatic technology in a Tecan Cavro product for the first time, this innovative module gives you an automated pipetting solution that is smaller than ever before, using air displacement to aspirate and dispense fluids. Correct operation is monitored by onboard pressure-based liquid level detection and self-diagnostics.

Excellent pipetting performance is demonstrated from 1-1000 µl, using a range of Tecan's high quality disposable tips, and an integrated tip ejector allows tips to be ejected by a simple command. The Cavro ADP's tip sensor can determine if a disposable tip has not been correctly picked up, or has fallen off the probe, flagging and logging errors to increase process security and aid validation of customer instrumentation, particularly for diagnostic applications.

To find out more about the Cavro Air Displacement Pipettor, visit www.tecan.com/components



Innovative HydroSpeed[™] plate washer launched

The HydroSpeed plate washer is an advanced system for optimized washing of cells, beads and ELISAs in 96- and 384-well formats. It offers full control over critical wash parameters, via an intuitive touchscreen interface, with extra gentle drop-wise dispensing and tunable aspiration settings to help avoid loss of material and maximize assay efficiency.

The HydroSpeed features advanced Cell Protection[™] settings for washing of adherent and loosely adherent cells, allowing the user to dial-in extra gentle wash parameters to suit their cells, microplates and application. The system's innovative Anti-Clogging[™] function takes the hassle out of ELISA washing by automatically rinsing and soaking the wash head when the instrument is idle between runs, and the Easy X-change system allows rapid removal and replacement of wash

heads for intense ultrasonic cleaning. The HydroSpeed uses a patent pending design with two magnets per well for high performance magnetic bead washing, offering fast bead settling and excellent recovery rates, and can also be equipped with a vacuum filtration module for processing of non-magnetic beads.

The HydroSpeed can be used in conjunction with Tecan's Connect[™] plate stacker for greater throughput efficiency, allowing semi-automated processing of up to 50 plates, or integrated into a Freedom EVO[®] liquid handling platform to offer fully automated processing for cell-, bead- and ELISA-based assay workflows.

To find out more on Tecan's HydroSpeed plate washer, visit www.tecan.com/cell-protection



HydroSpeed offers reproducible washing for magnetic bead-based assays



HydroSpeed plate washer

