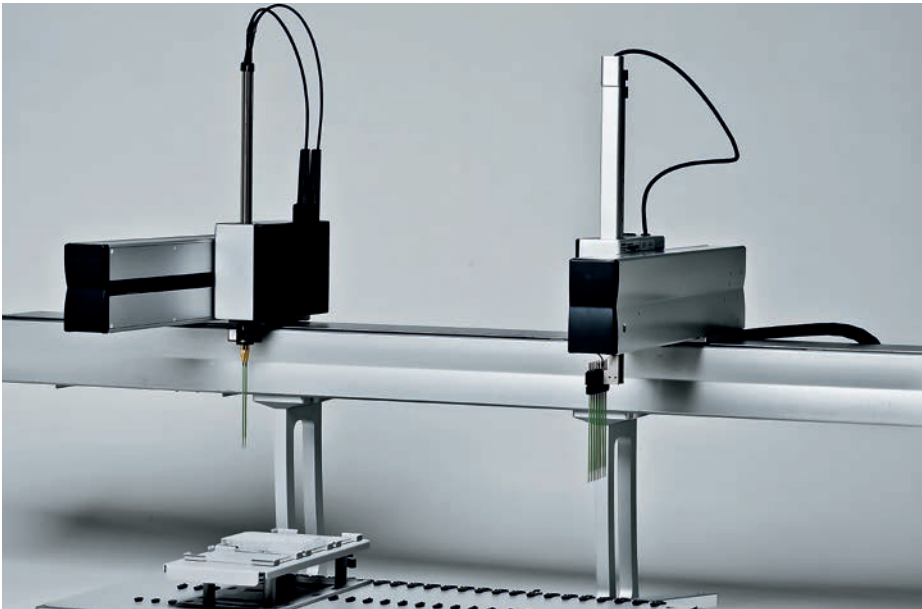
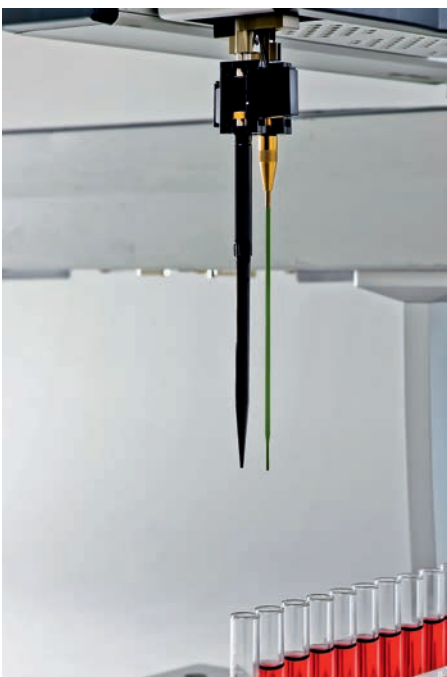


Extended flexibility of Cavro[®] Omni Robot



The Cavro Omni Robot offers precise OEM liquid handling

Tecan has introduced two extra options for the Cavro Omni Robot; the Dual Z and the ADP (Air Displacement Pipettor). Available with a broad choice of lengths, orientations and options for all three axes, the Cavro Omni Robot is designed to combine precision OEM liquid handling with straightforward integration. The launch



The Cavro Omni Robot Dual Z option can be fitted with a choice of fixed or disposable tips

of these additional pipetting modules provides instrument designers with even greater flexibility in instrument design.

The Cavro Omni Robot Dual Z option offers two independent pipetting channels in a single housing, doubling the robot's liquid handling capacity without having to mount an additional arm. This compact unit can be fitted with either fixed probes or disposable tip adapters, or a combination of both, with integrated capacitive liquid level detection (cLLD) and an adjustable pitch of 9 or 18 mm between channels. This elegant solution offers several potential advantages – especially where space is an issue – allowing instrument developers to increase throughput or perform additional operations without compromising on performance. For example, by fitting the Cavro Omni Robot Dual Z option with one fixed tip and one disposable tip adapter, it is possible to use the fixed probe for bulk reagent dispensing and disposable tips to pipette samples, avoiding the risk of cross-contamination while saving on consumables costs.

The Cavro Omni Robot is now also available with the Cavro ADP as an alternative to the existing liquid displacement pipetting technologies. Using pneumatic technology, this compact pipetting module has been

fully integrated and optimized for use with the Cavro Omni Robot's Universal Z arm, offering a number of additional benefits for OEM liquid handling. Its validated, maintenance-free design simplifies system integration, with on-board pressure-based liquid level detection (pLLD) and diagnostics to monitor correct operation. Together with the Cavro Omni Robot's standard cLLD, these functions provide exceptional pipetting performance and enhanced process security. As a self-contained, fully programmable pipetting option, the Cavro ADP is ideally suited to applications where access is restricted, and the module's integrated tip ejection option makes it possible for the Cavro Omni Robot to perform fully automated tip re-racking for the first time.

These additional pipetting modules further extend the versatility of the Cavro Omni Robot and, together with the system's extensive choice of options and finishes, provide exceptional flexibility for OEM liquid handling applications.

To find out more on Tecan's Cavro Omni Robot, visit www.tecan.com/omnirobot



The Cavro Omni Robot ADP option offers maintenance-free pipetting