

New Fluent™ features increase speed and flexibility

Spend more time doing the things that matter thanks to new Fluent laboratory automation solution* features designed to provide more streamlined and robust walkaway processing. These latest additions further extend the flexibility of Fluent, maximizing productivity and ensuring a precise match for your laboratory's automation needs.**

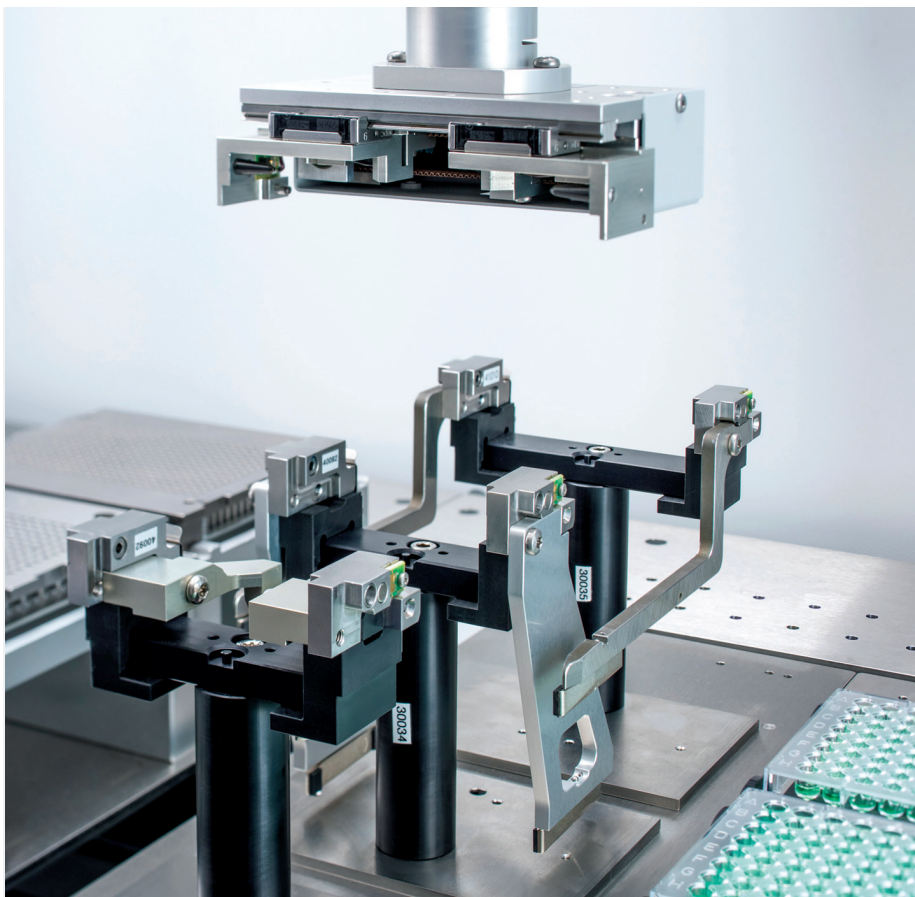
The new Finger Exchange System allows the workstation's Robotic Gripper Arm to automatically switch between gripper options – including Centric Fingers, Eccentric Fingers and Tube Fingers, plus customized options available through the Tecan Integration Group – on the fly. This ensures secure transport for virtually any automation-friendly plate or tube format, allowing straightforward sample reformatting and transfer of labware between centrifuges, storage modules,

processing devices and the workdeck, without the need for user intervention.

The flexibility of Fluent's high definition liquid handling has also been further enhanced, with the option to mount two Flexible Channel Arms on a single instrument. This pipetting option enables access to virtually any tube or well configuration, and provides a choice of air or liquid displacement pipetting technologies to suit your application.

Customers can now opt to combine both technologies on one platform, using air displacement for rapid pipetting with disposable tips, and liquid displacement for reagent distribution or bulk dispensing. This ensures you can always select the most appropriate pipetting technology for your application – offering the best of all worlds – and gives you the option to increase throughput by using both Flexible Channel Arms together to prepare a single plate or perform parallel analyses from common samples.

Understanding that confidence in workflow automation is critical to success, the Tecan team has also developed Method Recovery for FluentControl™. This powerful yet simple feature helps to minimize the disruption caused by an unexpected event – such as a power loss or computer failure – providing a straightforward, user-friendly way of salvaging a run. Using the integrated touchscreen, Method Recovery guides the operator through the steps required to re-establish an interrupted protocol, minimizing potential loss of valuable samples and reagents.



The new Finger Exchange System allows on-the-fly switching between gripper heads

To find out more about Tecan's Fluent workstations, visit www.tecan.com/fluent

* For research use only. Not for use in diagnostic procedures.

** New features due to be released soon.