

# Fluent™ – the new laboratory automation solution

Fluent is a unique automation concept built around the application-specific needs of laboratories, delivering more capacity and increased speed. Initially focusing on the complete automation of cell-based and biochemical assay workflows, Fluent brings together all the modules and devices required for automation of these assays in a single efficient and easy-to-use platform. It offers higher throughput, greater precision and improved reliability, allowing researchers to concentrate on what really matters – the results.

The Fluent solution for cell-based assays is designed to take the complexity out of cell biology research. Completely new from the ground up, it is available in three sizes – offering deck capacities from 30 to 72 microplates – to suit the throughput requirements of almost any laboratory. The patented Dynamic Deck™ uses a modular, multi-level design to offer exceptional deck capacity, allowing flexible integration of all the complementary devices required for cell-based research – including incubators,

washers and readers – into a single, fully automated system.

Developed to meet the everyday challenges faced by cell biology laboratories, Fluent aims to make cell-based assay workflows faster and more efficient. Liquid handling and labware logistics have never been easier, thanks to the instrument's three, task-specific arms which operate simultaneously to ensure timely completion of assays, minimizing the time cells spend outside of the incubator. Perfect coordination of the robotic arms is guaranteed by the platform's Path Finder™ move optimization technology, which continuously adapts to the changing topography of the workdeck, avoiding the need for laborious programming of individual arm movements, helping to accelerate both set-up and routine system use.

Even complex cell-based assays can be automated using the platform's FluentControl™ software, freeing researchers to focus on other tasks. Fluent's built-in touchscreen interface simplifies day-to-day

activities by guiding scientists through routine set-up and operation of the system for consistent, reproducible operation. It can provide graphical instructions and prompts for every step of the set-up process, from the number and type of cell plates required to the exact location of compound plates and assay reagents, helping to ensure compliance with SOPs and assay protocols. This is complemented by a PC-based programming environment which offers 'drag and drop' protocol development, 3D modeling and contextual checks to ensure rapid, easy scripting. FluentControl also offers straightforward integration and control of a wide range of Tecan and third-party devices, providing user-friendly, walkaway automation of virtually any cell-based assay workflow.

To find out more on Tecan's new Fluent laboratory automation solution, visit [www.tecan.com/fluent](http://www.tecan.com/fluent)

## Exceptional ease of use



Fluent's touchscreen interface guides users with clear graphical instructions and alerts for optimum simplicity and performance

## Unprecedented capacity, small footprint

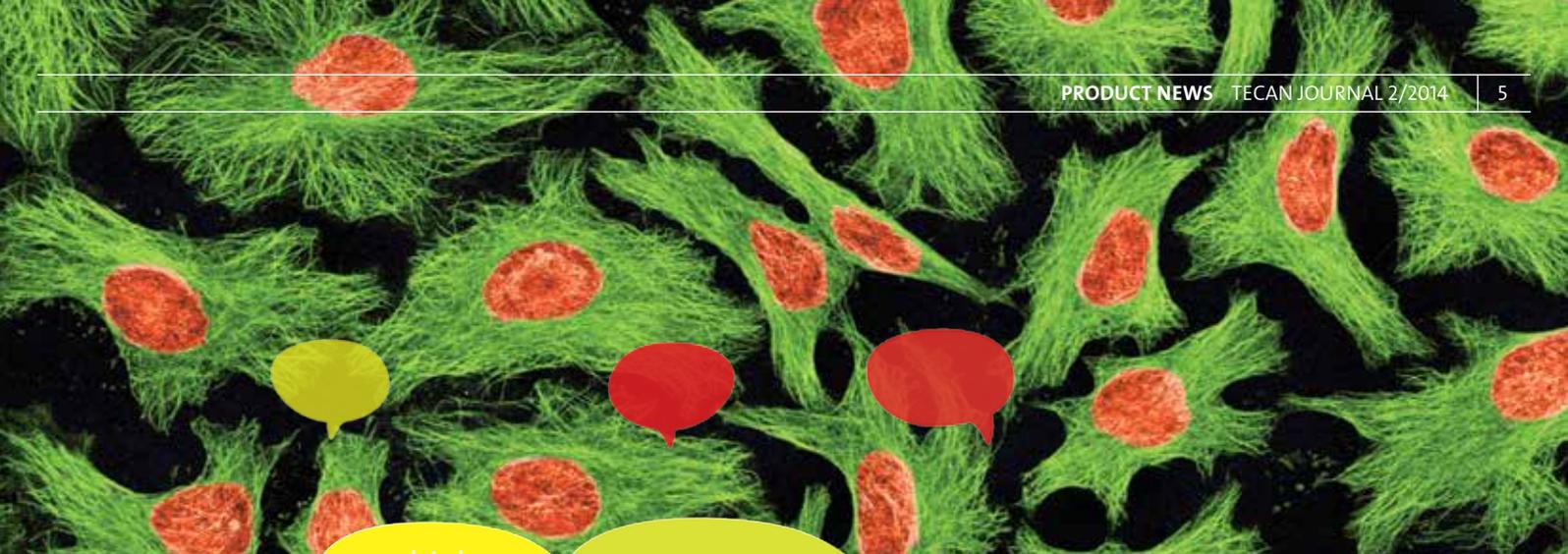


The Dynamic Deck offers unrivalled capacity and optimum efficiency for device integration

## Optimized cell protection



Simultaneous parallel movements of task-specific arms enhance speed and throughput



Improving productivity for complex cell-based research is a challenge. Processes are complicated, space is in short supply, lab personnel are confronted with more and more difficult tasks, and greater accuracy is always demanded. With the launch of the Fluent laboratory automation solution, all that changes. World-class productivity and simplicity are now within the grasp of every lab.

In the lab of the future, everybody will be fluent in cell-based assay automation

new

To see how Tecan is shaping the lab of the future, go to our webinar [www.tecan.com/webinars](http://www.tecan.com/webinars)

Watch the Fluent webinar now

