

# Added functionality with the MultiChannel Arm™ 384 gripper option

A gripper for the recently launched MultiChannel Arm™ 384 brings greater functionality to the high throughput pipetting head, and will be available as an option or upgrade. The cost-effective alternative to a robotic arm mounts onto the MultiChannel Arm 384, and shares its structure and x axis.

The gripper enables flexible access to all areas of the Freedom EVO® workstation with its unlimited 360-degree rotational axis, and the full range of its z axis allows the gripper to reach positions from the level of the worktable surface, up to

the highest shelf of a hotel. It is able to access objects located at the back of the instrument, such as labware stored in hotels and incubators, and can load/unload plates from readers or hotels on the right hand side of the platform. The gripper can maneuver a plate by grasping either its short or long side. The gripper also minimizes the risk of sample contamination by removing the plate lid only for the duration of a pipetting step, replacing the lid immediately afterwards.

*Scientific instrumentation. Not for use in human clinical or diagnostic procedures.*



## The Freedom EVOlyzer® takes routine clinical diagnostics tasks off your hands

Tecan has launched an enhanced version of the Freedom EVOlyzer® for fully guided, automated processing of ELISAs through its new generation Freedom EVOLution™ 2.0 software, which offers an exceptionally user-friendly graphical user interface with optional touchscreen and a number of new application-oriented features. The open-platform, IVD-D compliant Freedom EVOlyzer automates ELISA kits from a variety of manufacturers and manages the entire workflow, from organizing the initial worktable configuration to providing the final assay results.

The Freedom EVOLution 2.0 software performs many routine laboratory tasks, such as continuous sample loading and automated processing of pre-dilution factors. A volume pre-checking function ensures there are sufficient reagents for all scheduled tasks, allowing long walkaway times for the user with complete peace of mind. The software's

simple graphical loading dialogs and loading interface provide easy operation throughout the process, and its flexible scheduling options allow the order of assay processing to be defined and prioritized. Assays can also be bunched into profiles that facilitate everyday

routine tasks and the system is suitable for working in batch mode using worklists or via an ASTM connection.

For more information on Tecan's Freedom EVOlyzer, visit [www.tecan.com/evolyzer](http://www.tecan.com/evolyzer)

*The Freedom EVOlyzer*

