

# Three-in-one platform for automated filtration, separation and plate washing

Tecan's HydroFlex™, the compact 3-in-1 application platform for automated vacuum filtration, magnetic bead separation and microplate washing, provides increased productivity and reliable results by automating a range of 96-well format applications, including PCR clean-up and bead assays, ELISAs, cell assays and protein arrays. The HydroFlex automates key process steps such as

self-sealing of filtration plates, on-line vacuum control and fast dispensing of wash buffers, and offers a tunable vacuum range from -50 mbar to -850 mbar for both high and low vacuum applications.

The HydroFlex is easily operated as a stand-alone instrument with its built-in keypad, or by using the intuitive

HydroControl™ software, which has been designed to meet the FDA CFR 21 part 11 regulations. The platform can be integrated with Tecan's Freedom EVO® liquid handling workstations or Freedom EVOlyzer® ELISA analyzer.



*The HydroFlex platform for automated vacuum filtration, magnetic bead separation and microplate washing is available in various colours*