

Automation doesn't get easier than this!

Tecan's Smart Automation solutions offer straightforward liquid handling for new users

Tecan is launching a range of benchtop solutions designed to provide fast, simple and cost-effective automation of liquid handling tasks. Based on the Freedom EVO® 75 platform, these preconfigured systems are designed to eliminate tedious manual liquid handling tasks without the need for previous automation experience.

Each Smart Automation solution is supplied with optimized, ready-to-run protocols, and includes all of the modules and ancillary devices necessary to automate your manual processes.

The first Smart Automation solutions to be launched are:

Pickolo™ colony picking; for fully automated colony picking

NucleoSpin® purification; for fully automated, vacuum-based nucleic acid extraction using MACHEREY-NAGEL NucleoSpin kits

NucleoMag® purification; for fully automated, magnetic bead-based nucleic acid separation using MACHEREY-NAGEL NucleoMag kits.



A dedicated user interface and easy-to-follow application guide ensure that even inexperienced users can get their systems up and running quickly, with touchscreen operation also available for some applications. These complete solutions provide an ideal introduction to the throughput, precision and reproducibility offered by Tecan's liquid handling systems, with the flexibility to expand and adapt as your laboratory needs change.

To find out more on Tecan's Smart Automation solutions, visit www.tecan.com/smartautomation

Tecan and Promega offer real-time cytotoxicity analysis

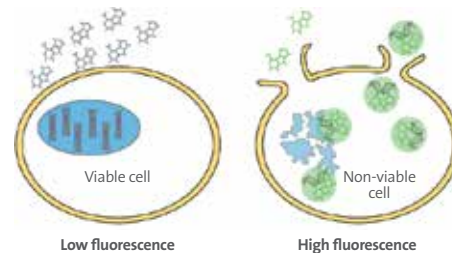


Tecan has worked with Promega to demonstrate the suitability of the Infinite® 200 PRO and HP D300 Digital Dispenser for real-time kinetic investigation of cytotoxicity using Promega's CellTox™ Green Dye.

The CellTox Green Cytotoxicity Assay measures changes in membrane integrity that occur as a result of cell death. This non-toxic dye allows real-time fluorescence-based measurements for up to 72 hours, and is ideally suited to mode-of-action studies. Combining the CellTox Green assay with the advanced environmental control and cell biology

oriented functions of the Infinite 200 PRO – including the Gas Control Module (GCM™), Optimal Read (OR), linear and orbital shaking, temperature control and automated z-focusing – offers an unparalleled solution for long-term kinetic studies of cytotoxicity. By providing stable culture conditions within the reader chamber, this set-up allows continuous measurement over the duration of the experiment, ensuring high quality data without the need for manual intervention or transfer of microplates between an incubator and the reader.

The assay workflow can be further improved by use of the HP D300 Digital Dispenser to provide on-demand assay plate set-up. Offering non-contact dispensing from 13 pl to 10 µl, this innovative instrument allows small molecules in DMSO to be dispensed directly into the assay plate, eliminating the need for time-consuming and wasteful serial dilutions. With a range of features specifically tailored towards cell-based and drug discovery applications – including plate shaking to



The CellTox Green Dye is excluded from viable cells, but binds to DNA from cells with compromised membrane integrity, resulting in an increase in fluorescence

prevent a bolus of DMSO from sinking onto plated cells, automatic plate randomization and straightforward generation of drug-drug interaction studies – the HP D300 is designed to increase productivity with minimal set-up time for cell-based and biochemical studies.

To find out more on Tecan's Infinite 200 PRO and GCM, visit www.tecan.com/gcm

To learn more about Direct Digital Dispensing, go to www.tecan.com/digitaltitration