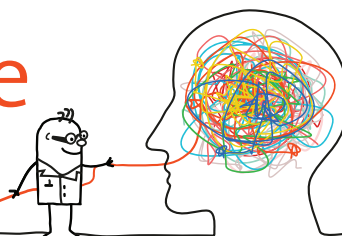


Software updates enhance flexibility of HP D300



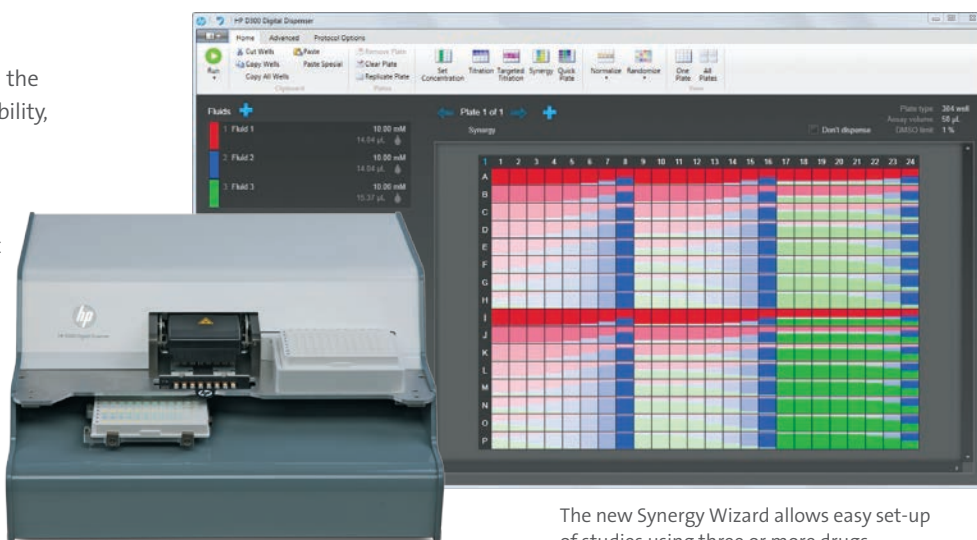
The latest software updates for the HP D300 Digital Dispenser ensure even greater versatility for drug discovery applications. The recently released version 2.2 software includes a number of updates to improve day-to-day operation of the HP D300 for new and existing customers, making it even faster to set up and run your experiments.

Format flexibility has been increased with the addition of 1,536-well microplate compatibility, complementing the instrument's existing 12- to 384-well plate direct dispensing capabilities. This high density format is particularly suited to assay development applications, providing better integration with downstream activities by allowing experiments to be designed and optimized in the same labware type used throughout the workflow. The instrument's Synergy Wizard has also been improved, allowing straightforward investigation of drug-drug interactions

between multiple compounds of interest. Ideal for cutting-edge research, the new Synergy Wizard allows the user to quickly and easily set up synergy studies for three or more drugs, developing complex experiments with ease. Alongside a host of minor updates to simplify instrument use, these

software upgrades are designed to accelerate experimental set-up for a wide range of drug discovery activities.

To find out more about the HP D300 Digital Dispenser, visit www.tecan.com/digitaltitration



The new Synergy Wizard allows easy set-up of studies using three or more drugs

Tecan Academy to be available worldwide

Tecan's online training tool, the Tecan Academy, will be launched globally in 2014, offering basic training and certification for new operators. The Academy's library of online content is continuously updated with new material, with recent additions including a new certification course for Freedom EVO 75 operators and a 'How to' guide on using the Infinite® series of multimode readers for fluorescence measurements. Each distinct learning module is designed to enhance user understanding for reduced operational errors and higher productivity, all without having to leave your desk!

Many users throughout Europe and the US are already enjoying the benefits of the greater instrument knowledge offered by the Tecan Academy. Ben Davis, a Tecan

Academy user and Scientist in the Technical Development department at LGC Genomics, commented: "The Tecan Academy was recommended to us as a user-friendly way of improving day-to-day operation of our Tecan multimode reader for QC testing. It is very easy to use and navigate, with a clean design and concise graphical representations of the technical aspects of the instrument. Using the Academy has certainly provided a useful refresher on how to apply theoretical understanding to practical investigations, and it was very convenient to use, allowing me to log in and out as required."

To find out more on the Tecan Academy, visit www.tecan.com/academy



The Tecan Academy offers convenient, web-based training