

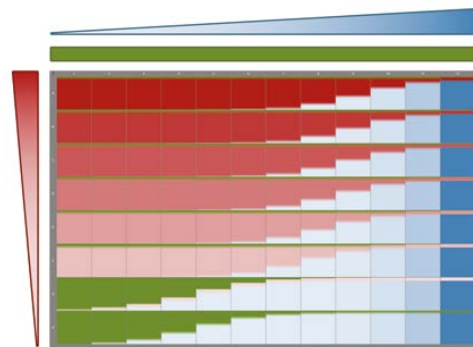
HP D300 now offers aqueous liquid dispensing

The popular HP D300 Digital Dispenser now allows customers to titrate low volumes of biomolecules in aqueous solutions, offering greater flexibility for drug discovery workflows. Users can choose between dispensing DMSO, for small molecule studies, and aqueous solutions in combination with a surfactant, for the investigation of proteins, antibodies, enzymes and nucleic acids.

Newly launched T8+ and D4+ Dispenseheads, combined with updated software, allow both new and existing customers to benefit from this breakthrough, offering straightforward set-up of dose-response curves and synergy studies. New 'fluid class' options allow users

to optimize liquid dispensing performance to their specific buffers and surfactants, and an Enzyme Profile function enables fast and easy set-up of titrations using differing combinations of enzymes, substrates and other compounds. These innovative features, combined with the system's proven low volume dispensing performance, make it easy to perform complex experiments – such as enzyme characterization, K_M determination and exploring the effects of enzyme inhibitors – ensuring a comprehensive solution for drug discovery applications.

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



■ Enzyme - Fixed concentration
■ Substrate - Titrated
■ Compound - Titrated

The new Enzyme Profile function enables rapid set-up of any combination of enzymes, substrates and compounds with just a few clicks

Cavro® Disposable Tips launched for OEM customers

Tecan has created a range of Cavro Disposable Tips for use with the Company's popular Cavro brand OEM components. Offering customers a selection of premium consumables to complement their Cavro Omni Robot and Cavro Air Displacement Pipettor (ADP) options, these high quality tips have been developed and validated to ensure optimal performance and reliable operation.

All Cavro Disposable Tips are manufactured to the same high quality standards as Tecan's broad portfolio of consumables for life sciences customers, using state-of-the-art production processes and integrated quality control procedures to ensure consistent pipetting performance and provide complete peace of mind. The conductive polypropylene tips are available in a variety of tip volumes, and can be supplied in customer-branded packaging for large volume orders, providing a complete liquid handling solution to enhance your instrument design.

To find out more on Tecan's Cavro Disposable Tips, visit www.tecan.com/cavrodit



Cavro Disposable Tips offer exceptional performance and peace of mind