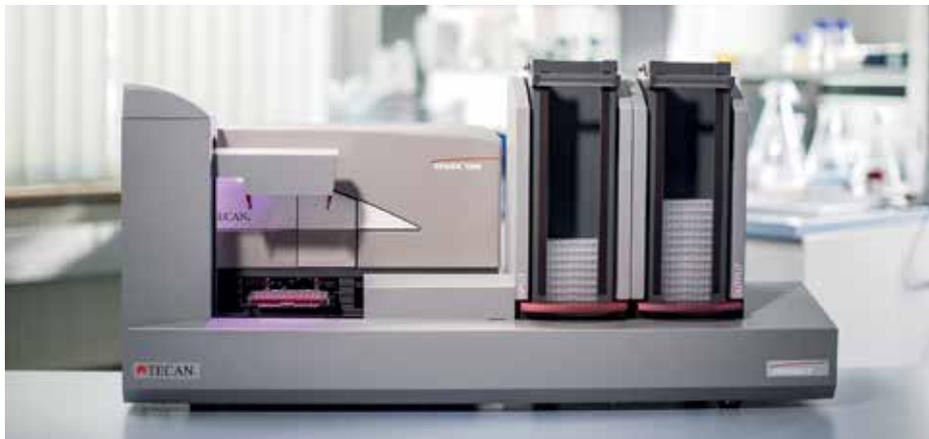


Walkaway batch processing for the Spark™ 10M

The Spark 10M multimode reader can now be combined with the Connect™ microplate stacker to offer semi-automated batch processing of up to 50 assay plates. A newly developed software interface allows seamless operation of the Connect stacker through the reader's SparkControl™ software, helping to streamline your laboratory workflow for greater productivity.

The Connect is a versatile microplate stacker designed to offer fast, smooth transfer of plates to and from the host instrument. Already popular with customers using Infinite® series readers and HydroSpeed™ plate washers, this proven and reliable system can now be used in conjunction with the powerful Spark 10M reader. All common microplates conforming to ANSI/SLAS standards – from 6- to 1,536-wells – can be used with Connect, with a choice of



The Connect stacker offers walkaway processing for up to 50 microplates using the Spark 10M

30- or 50-plate cassettes to suit a variety of throughputs. Optional dark covers are also available to protect light sensitive reactions, making the system ideally suited to a wide range of applications in genomics, proteomics and drug discovery.

To find out more about Tecan's Connect stacker and Spark 10M multimode reader, visit www.tecan.com/connect and www.tecan.com/spark

Tecan's D300e offers fast liquid transfer down to picoliter levels

Tecan has introduced the D300e Digital Dispenser* for applications that need faster, reliable dispensing down to picoliter volumes. This innovative, simple-to-use device is ideal for the creation of assay plates, and can cut set up times from hours or days to just minutes. Users can choose between dispensing aqueous solutions in combination with a surfactant – for the investigation of proteins, antibodies, enzymes and nucleic acids – and DMSO, for small molecule studies, offering greater flexibility for life sciences workflows.

The system's sophisticated yet easy-to-use software guides you through every operation, minimizing training times and problem-solving call-outs, and providing straightforward set up of enzyme profiles, dose-response curves and synergy studies. Even the most complex experimental plate layouts – such as drug-drug interaction studies – can be quickly and easily generated, with optional randomization to

reduce the impact of edge effects and increase data integrity. The D300e is compatible with a wide range of microplate formats from 12 to 1,536 wells, including deep-well plates and, for maximum flexibility, can even accommodate ANSI/SLAS-format tube racks.

To find out more about Tecan's D300e Digital Dispenser, visit www.tecan.com/D300e

*For research use only. Not for use in diagnostic procedures.



Tecan's D300e Digital Dispenser