

Tecan and *p*ION offer new automated drug screening solution



*p*ION's Double-Sink PAMPA integrated with Tecan's Freedom EVO platform

Following on from the successful launch of manual assay systems for permeability and solubility, Tecan and *p*ION INC have co-developed an automated solution for drug permeability and solubility screening. This new solution is both scalable and upgradable according to the user's changing requirements, from the entry level manual technique to a fully automated system for high throughput applications.

*p*ION's Double-Sink™ parallel artificial membrane permeability assay (PAMPA) helps to identify the most promising lead compounds well in advance of cell-based methods, and has been integrated with the Freedom EVO® workstation and Infinite® M200 plate reader in the PAMPA Evolution™ system. This platform offers enhanced

walkaway time and high throughput for fully automated processing of permeability assays, while automated miniaturized shake-flask solubility assays can be done with the similarly configured μ SOL Evolution™. The flexibility of the Freedom EVO workstation also allows a single instrument to perform both permeability and solubility assays, as well as other ADME assays, in a combined set-up. The advanced Quad4 Monochromators™ technology in the Infinite M200 provides excellent sensitivity, and an optional MultiChannel Arm™ 96 further increases throughput with rapid sample transfers.

“Over the years, we have invested heavily in developing new applications for our patented Double-Sink PAMPA, allowing our

licensed customers to conduct cost-effective screening prior to the more expensive cellular and *in vivo* GIT and BBB testing,” explained Dr Alex Avdeef, *p*ION CSO and President. “We are delighted to continue working with Tecan, with its Freedom EVO platform demonstrating once again the consistently high quality and reliability of all its products.”

To find out more on Tecan's PAMPA Evolution system, visit www.tecan.com/pampa

Double-Sink, PAMPA Evolution and μ SOL Evolution are trademarks of *p*ION INC, Woburn, USA.