

# Direct Digital Titration triumphs at SLAS2012

The first annual SLAS Conference and Exhibition held in San Diego, California, in February heralded success for Tecan, with the HP D300 Digital Dispenser receiving the SLAS2012 New Product Award Designation.



Jessica Merlino, Tecan Product Manager, and Christie Dudenhoefer, HP Instrument Design Lead, accept the award.



An intensive selection process involved an initial assessment phase, where individual judges visited their allocated booths to evaluate and rank each entry before coming together to establish a short list of candidates for the award. Phase two saw the team of judges rating the top nine entries against specific, detailed criteria and, of more than 60 new products launched at SLAS2012, the innovative HP D300 was one of just four chosen to receive this prestigious award.

And it wasn't just the judges who were impressed by the HP D300. The instrument received widespread interest from delegates, who were impressed by its user-friendly interface and ability to rapidly deliver picoliter to microliter volumes of drug compounds. This latest award is further proof of the convenience and flexibility of Direct Digital Titration and its potential to improve productivity and accelerate the discovery of better drugs.

To find out more about Direct Digital Titration, go to [www.tecan.com/digitaltitration](http://www.tecan.com/digitaltitration)

## 2012 Tecan Symposium – mass spectrometry

Boston, October 2012

Tecan is proud to announce its fifth Tecan Symposium, to be held this year from the 23<sup>rd</sup> to the 25<sup>th</sup> of October in Boston, USA. This year's topic will be *'Mass spectrometry – the expanding role in life sciences and diagnostics'* and will include: the application of mass spectrometry (MS) in diagnostics and other fields; the role of mass spectrometry in basic research; the collection, preparation and storage of samples for MS; and novel and emerging MS technologies. As in previous years, attendance at the symposium is limited to a small cohort of scientists in order to encourage plenty of networking opportunities among thought leaders in the area of mass spectrometry.

For more information, visit [www.tecan.com/symposium](http://www.tecan.com/symposium)

