

Automated mass spectrometry sample preparation; the key to an efficient workflow

In recent years, monumental advances have been made in mass spectrometry (MS) instrumentation, but sample preparation has not enjoyed the same rate of development, making it a major bottleneck in the laboratory workflow. The issues associated with manual processing have hindered the uptake of MS innovation in the life science industry. To help overcome these limitations, Tecan offers Freedom EVO-based end-to-end process automation solutions for even the most challenging sample preparation protocols, liberating scientists from the tedium of manual sample preparation.

A wide range of protocols – ranging from the low complexity 'dilute and shoot' and protein crash methods used in toxicology workflows, to the more complex solid phase (SPE), liquid-liquid (LLE) extraction or protein purification/digestion techniques – are routinely used for MS sample preparation. Automation of these processes is key to meeting the high throughput and quality demands of customers. Tecan's freely-configurable, automated extraction and purification platforms eliminate manual errors, allowing staff to focus on downstream MS analysis, streamlining laboratory workflows

and enhancing productivity. For common MS sample preparation applications, the easy-to-use TouchTools™ interface places solutions at customers' fingertips, making automation easier than ever before.

To discover how automated MS sample preparation is benefitting ADPEN Laboratories, see pages 16-17.

To find out more on Tecan's MS sample preparation solutions, visit www.tecan.com/lcms

Go configure! Receive a customized quote fast!

Customers can now plan and create over 100 different designs of Cavro® Omni Robot using Tecan's new online graphical Cavro Omni Robot Configurator. This straightforward, web-based application – available for PCs or iPads® – provides laboratory instrument designers with a quick and easy way to explore the various options available for the Cavro Omni Robot, helping to accelerate their product development.

The Cavro Omni Robot Configurator details the complete range of standard options for the Cavro Omni Robot, including the various axis sizes and arrangements, arm functions and liquid handling configurations, as well as a number of finishing features, such as covers and end caps. Users can generate an interactive 3D model of a selected Cavro Omni Robot specification that updates immediately as different options are chosen, and the image can be rotated on the screen to help visualize the finished robot from different perspectives. This provides engineers with a visual confirmation of how the selected

configuration will fit, as well as quick and easy verification of the dimensions that might affect instrument design. Once the required specification has been selected, the Cavro Omni Robot Configurator generates a complete parts list which can be printed or e-mailed directly to the Tecan sales team for a quote.

To build your Cavro Omni Robot now, go to www.tecan.com/cavroomniconfigurator

