

# Freedom EVOLyzer<sup>®</sup> received with great enthusiasm in Spain



The Freedom EVOLyzer, Tecan's new generation of automated ELISA systems, has proved to be increasingly popular with laboratories throughout Spain since being introduced to the country in July 2006 by Inverness Medical Iberica (IMI), one of Tecan's principle diagnostic distributors for the area. IMI, formerly part of Innogenetics Diagnostics, has a long and proven track record as a distributor for professional diagnostic products in Spain, with a strong focus on enzyme-linked immunosorbent assays (ELISAs) for infectious diseases, diagnostics and autoimmunity. IMI's customers range from large hospital laboratories performing routine testing, to reference laboratories and testing laboratories for small- to medium-sized hospitals.

Tecan's cooperation with IMI started in the early 1990s and, during the last decade, IMI has placed a broad

range of Tecan instruments in Spain, including Genesis RMP™ systems, Minilyzers, Profiplot™ analyzers, Sunrise™ plate readers and Columbus™ plate washers. IMI is a distributor for nine international reagent manufacturers and the company's combined kit portfolio consists of more than 300 different tests, which IMI has successfully automated using Tecan's open ELISA platforms such as the Minilyser and the Genesis RMP. The addition of the Freedom EVOLyzer has significantly strengthened IMI's IVD-D compliant portfolio.

"We chose to take full advantage of the large range of customization and configuration possibilities that the Freedom EVOLyzer offers, such as being able to freely combine stainless steel tips and disposable tips on the same instrument," explained Mr Sergio Gordillo, Automation Product Manager at IMI. "The instrument's flexibility for both hardware and software options means

we are able to satisfy almost all kinds of ELISA customer with just one single platform, which has been a key factor in the Freedom EVOLyzer's success. We always deliver these instruments with a touchscreen monitor, because the Freedom EVOLyzer software has a touchscreen compatible interface available in both Spanish and Catalan, and this has proven to be a real hit among our local customers."

"We have been able to transfer 100% of our kits to the Freedom EVOLyzer with excellent results," continued Mr Gordillo. "This new system is a more than worthy successor of Tecan's Genesis RMP and its function scope shows a clear improvement in terms of both throughput and flexibility."

"We have relied on Tecan instrumentation for a good number of years now and it has proven to be the right choice. With the Tecan instruments and our broad range of assays we have managed to gain a significant market share of the Spanish market and, with more than 80 ELISA platforms placed in Spain, everybody knows that we are a force to be reckoned with!" claimed Mr Gordillo. "With the Freedom EVOLyzer we can now offer superior kits, automated with superior instruments supported by superior service. What customer would ever say no to that?"

*The Freedom EVOLyzer for automated ELISAs*

