

# Automated multiplex assays offer novel newborn screening solution

## Luminex®

Tecan is collaborating with Luminex® Corporation to develop an innovative, automated newborn screening solution designed to rapidly identify potential health issues in infants. Luminex's Luminex 200™ instrument with xMAP® technology is being integrated with Tecan's Freedom EVO® liquid handling platform and HydroFlex™ washer to create the 996x™.

xMAP technology allows large numbers of biological tests to be performed and analyzed simultaneously, and xMAP newborn assays offer high throughput testing for multiple biomarkers linked to critical diseases. Screening enables early detection of potential health problems, allowing effective treatment to be initiated, and the 996x will provide a walkaway automated solution for cost-effective, multiplexed newborn screening. This new platform, which is expected to be launched in 2011, will help laboratories enhance efficiency, save vital resources and analyze more samples, more quickly.

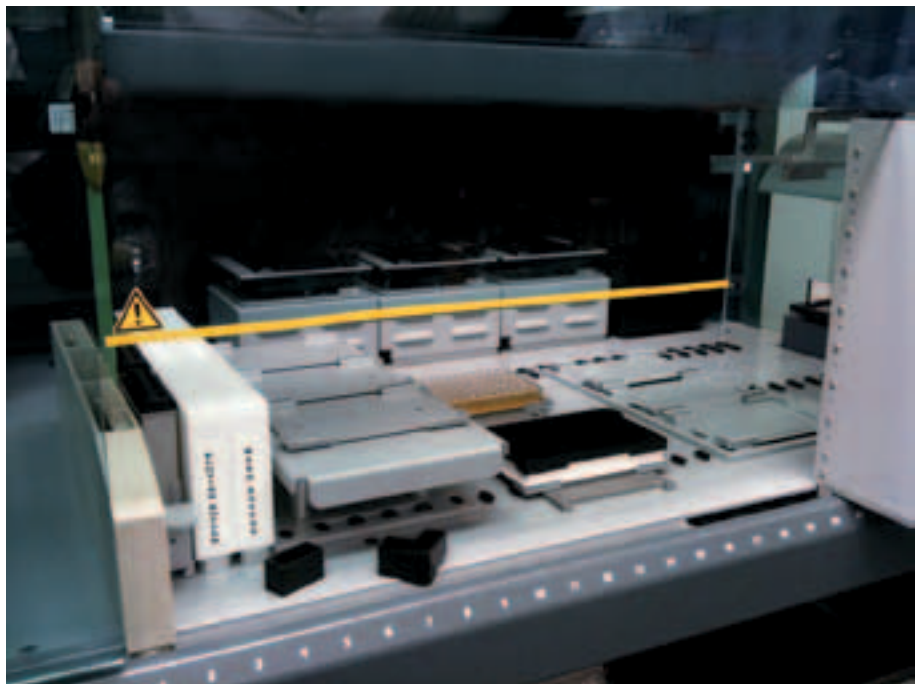
To find out more about Tecan's Freedom EVO workstations, visit [www.tecan.com/freedomevo](http://www.tecan.com/freedomevo)

To find out more on Luminex Corporation or xMAP technology, visit [www.luminexcorp.com](http://www.luminexcorp.com)

Luminex 200 and 996x are trademarks, and Luminex and xMAP are registered trademarks of Luminex Corporation.



Project team members from Luminex and Tecan (left to right): Kyle Knight (Luminex), Renate Kehrli and Christoph Beck (Tecan)



xMAP technology integrated onto Freedom EVO platform to create 996x