

Tecan Group and Luminex Corporation Partner to Develop Automated Newborn Screening Solution

New solution combines multiplexing and automation technologies to create novel system that provides faster results and saves valuable time and resources in newborn screening laboratories

Männedorf, Switzerland, and Austin, Texas – October 12, 2010 – The Tecan Group (SIX Swiss Exchange: TECN) and Luminex Corporation (NASDAQ: LMNX) have announced that they will collaborate on the development of a new automated newborn screening solution that will provide faster results for infants' potential health concerns. The companies have entered into a global OEM agreement that allows them to combine Luminex® xMAP® technology and LX 200™ instrument with Tecan's flexible and reliable Freedom EVO® liquid handling platform and HydroFlex™ washer. The new automated newborn screening solution will help laboratories enhance efficiency, save vital resources and analyze more samples more quickly. Luminex expects to launch the automated newborn screening solution in 2011.

Thomas Bachmann, CEO of Tecan, said: "We are delighted that we have been chosen by Luminex to supply a fully automated solution for such an important application. This global OEM agreement is building on a long, well established partnership of creating the most advanced automated solutions in the research market for Luminex xMAP technology."

Luminex xMAP technology enables large numbers of biological tests to be conducted simultaneously and analyzed quickly, cost-effectively and accurately. Once it receives regulatory clearance, NeoPlex4™ will be a fully automated, walk-away testing platform for multiplexed newborn screening. These high-volume screening assays test newborn infants for several different biomarkers linked to critical diseases. Each of these conditions can cause severe mental and physical disability if untreated. If detected early through testing, effective treatment can be initiated. Multiplexing these tests together from one sample offers significant improvements in testing efficiency.

Patrick Balthrop, CEO of Luminex, said: "Automation continues to be a top priority for Luminex and our customers. The high testing volumes in the newborn screening market require full sample-to-answer automation for the thousands of specimens processed by each customer per day. We evaluated several different robotic solutions to achieve this full automation for our xMAP technology and determined that Tecan was the best option available for this application. We have had a very effective relationship with Tecan and we are proud to be partnered with them in this effort."



2

About Tecan

Tecan (www.tecan.com) is a leading global provider of laboratory instruments and solutions in biopharmaceuticals, forensics, and clinical diagnostics. The company specializes in the development, production and distribution of instruments and automated workflow solutions for laboratories in the life sciences sector. Its clients include pharmaceutical and biotechnology companies, university research departments, forensic and diagnostic laboratories. As an original equipment manufacturer, Tecan is also a leader in developing and manufacturing OEM instruments and components that are then distributed by partner companies. Founded in Switzerland in 1980, the company has manufacturing, research and development sites in both Europe and North America and maintains a sales and service network in 52 countries. In 2009, Tecan generated sales of CHF 392 million (USD 361.2 million; EUR 259.6 million). Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TK: TECN/Reuters: TECZn.S/ ISIN CH0012100191).

About Luminex Corporation

Luminex Corporation develops, manufactures and markets proprietary biological testing technologies with applications throughout the diagnostic and life sciences industries. The Company's xMAP[®] multiplex solutions include an open-architecture, multi-analyte technology platform that delivers fast, accurate and cost-effective bioassay results to markets as diverse as pharmaceutical drug discovery, clinical diagnostics and biomedical research, including the genomics and proteomics markets. The Company's xMAP Technology is sold worldwide and is already in use in leading clinical laboratories as well as major pharmaceutical, diagnostic and biotechnology companies. Further information on Luminex Corporation or xMAP Technology can be obtained at www.luminexcorp.com.

For further information:

Tecan Group

Martin Braendle
Head of Corporate Communications & Investor Relations
Tel. +41 (0) 44 922 84 30
Fax +41 (0) 44 922 88 89
investor@tecan.com
www.tecan.com

Luminex Contacts:

Luminex Corporate Contact:
Harriss T. Currie, Vice President, Finance and Chief Financial Officer
512.219.8020
hcurrie@luminexcorp.com

Matthew Scalo
Sr. Director Investor Relations
512.219.8020
mscal@luminexcorp.com

Luminex Media Contact:
Aaron DeLucia
512.241.2249
aaron.delucia@porternovelli.com