

## **Change in the Tecan Board of Directors**

- **Long-standing Board of Directors Chairman Mike Baronian will not stand for re-election at the ordinary Annual General Meeting 2009**
- **The Board will propose the current members Heinrich Fischer, Gérard Vaillant, Prof. Dr. Peter Ryser and Dr. Juerg Meier for re-election**
- **Rolf A. Classon, Dr. Lukas Braunschweiler and Dominique F. Baly will be proposed for election as new members**
- **Rolf A. Classon expected to take over as Chairman**

**Männedorf, Switzerland, March 4, 2009** – Change in the Board of Directors of Tecan (SIX Swiss Exchange: TECN): in line with a periodic renewal, the Tecan Board of Directors proposes to the ordinary Annual General Meeting on 22 April 2009 Heinrich Fischer, Gérard Vaillant, Prof. Dr. Peter Ryser und Dr. Jürg Meier for re-election and Rolf A. Classon, Dr. Lukas Braunschweiler and Dominique F. Baly as new members of the Board. Rolf Classon is expected to succeed Mike Baronian and take over as Chairman.

In the name of the Board of Directors, Heinrich Fischer, Vice Chairman, thanked the Board members who will resign at the Annual General Meeting Mike Baronian, Cleto de Pedrini and Armin Seiler “for their long-standing loyal and successful contribution. They played an important role for Tecan which is now well positioned for the future.”

### **Three new members will be proposed**

Rolf A. Classon, born 1945, is on the Board of various US and European companies in the life science sector. Up until 2006 he held various top management positions at large international corporations in the pharmaceutical sector, including CEO at Bayer Healthcare and CEO of Hillenbrand Industries. A Swedish citizen, he holds a degree in chemical engineering from the University of Gothenburg.

Dr. Lukas Braunschweiler, born 1956 and Swiss citizen, graduated from the Swiss Federal Institute of Technology (ETH Zurich) in physical chemistry. He was appointed CEO of Ruag Holding AG in December 2008 and will take up this role in June 2009. From 2002 to date he has been Delegate of the Board of Directors and CEO of Dionex, a listed US-based

manufacturer of laboratory instruments. He previously held executive positions at a number of major high-tech companies including Mettler-Toledo and Landis & Gyr.

Dominique F. Baly, born 1949, graduated from the University Louis Pasteur in Strasbourg, France, in Pharmacy. He is a French and US citizen. From 1972 until 2008, Dominique F. Baly held various top management positions at Millipore Corporation in Europe and in the US, including President of the Bioscience Division, President of the Laboratory Water Division and Vice President & General Manager, Millipore Europe.

### **About Tecan**

Tecan ([www.tecan.com](http://www.tecan.com)) is a leading global provider of laboratory instruments and solutions for the biopharma, forensic and diagnostic industries. The company specializes in the development, production and distribution of automation solutions for laboratories in the life sciences sector. With its subsidiary REMP ([www.remp.com](http://www.remp.com)), Tecan is the market leader in automated laboratory storage and logistics systems. Its clients include pharmaceutical and biotechnology companies, university research departments and diagnostic laboratories. Founded in Switzerland in 1980, the company has manufacturing, research and development sites in both North America and Europe and maintains a sales and service network in 52 countries. Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TK: TECN/Reuters: TECZn.S/Swiss security number: 1210019).

### **For further information please contact:**

#### **Tecan Group**

Mike Baronian

Chairman of the Board of Directors

+41 31 810 40 10

[www.tecan.com](http://www.tecan.com)

