

## **Tecan Names New Head of Business Unit Sample Management**

- **Domingo Messerli named Head of Business Unit Sample Management (formerly REMP) and a member of the Executive Board**
- **A manager with international experience in the industry**

**Männedorf, Switzerland, June 23, 2008** – The Board of Directors of Tecan Group AG (SWX: TECN) appointed Domingo Messerli Head of Business Unit Sample Management (formerly REMP) and a member of the Executive Board. He will assume the position at Tecan on July 1<sup>st</sup>, 2008. Since the end of March 2008 the Business Unit was temporarily managed by Thomas Bachmann, CEO of the Tecan Group. With its Business Unit Sample Management Tecan is the premier supplier of automated laboratory storage and logistics systems.

From 2005 until spring 2008, Domingo Messerli was CEO of Elektrizitäts AG in Basel, Switzerland, a provider of electrical solutions and services. Before that he held a variety of management positions at Glatt AG in Pratteln, Switzerland, an international manufacturer and distributor of complex capital goods for the pharmaceutical, food and fine chemical industries. For four and a half years Domingo Messerli was the CEO of the Swiss headquarters of the Glatt Group as well as its associated international subsidiaries. In addition he was the Head of Global Sales and Marketing for the company's Business Unit Pharmaceuticals. Domingo Messerli holds a degree in Mechanical Engineering and an Executive MBA from the University of St. Gallen. He is 49 years old and Swiss citizen.

Thomas Bachmann, CEO of Tecan, comments as follows: "We are pleased to have Domingo Messerli take over at the helm of our Sample Management business. He is a manager with international experience and a broad range of expertise in the industry. His job at Tecan will be to make efficient use of the potential for growth in Sample Management and improve the unsatisfactory profitability of that business unit. We are confident that Domingo Messerli will work energetically to advance our Sample Management business and will make a valuable contribution to the prosperity of the Tecan Group."

### **About Tecan**

Tecan ([www.tecan.com](http://www.tecan.com)) is a leading global supplier 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 life science laboratories. Through its REMP subsidiary ([www.remp.com](http://www.remp.com)), Tecan is the premier supplier of automated laboratory storage and logistics systems. Tecan's clients include pharmaceutical and biotech companies, university research



2

departments, diagnostic laboratories and companies as well as forensics laboratories. Founded in Switzerland in 1980, the company has production, research and development sites in both North America and Europe and maintains a sales and service network in 52 countries. In 2007, Tecan achieved sales of CHF 414.4 million (USD 345.3 million; EUR 252.7 million). Registered shares of the Tecan Group are traded on the SWX Swiss Exchange (TK: TECN/Reuters: TECZn.S/Valor: 1210019).

**For further information, please contact:**

**Tecan Group**

Dr. Rudolf Eugster  
Chief Financial Officer

[investor@tecan.com](mailto:investor@tecan.com)  
[www.tecan.com](http://www.tecan.com)

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