

## Tecan Announces Change in Management Team

- **Matthias DÜchting, regional head of sales and member of the Group Executive Committee is leaving Tecan**

**Männedorf, Switzerland, November 18, 2009** – The Tecan Group (SIX Swiss Exchange: TECN) today announced a change in its Management Team. Matthias DÜchting, Head of the European sales organization and member of the Group Executive Committee, will leave Tecan by mutual agreement. Jürg Dübendorfer, Head of Business Unit Services & Consumables, takes over the responsibility for the Sales Region Europe on an interim basis with immediate effect until a successor is found. Jürg Dübendorfer has been with Tecan since 2001 and has been a member of the Group Executive Committee since 2006.

Matthias DÜchting joined Tecan in October 2008 and was subsequently appointed as Head of sales for the "Europe & NEMEA" region, which covers the markets in Europe, Russia, the Middle East, Africa and India.

### 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. In 2008, Tecan generated sales of CHF 396 million (USD 366.7 million; EUR 250.7 million). Registered shares of Tecan Group are traded on the SWX Swiss Exchange (TK: TECN/Reuters: TECZn.S/Swiss security number: 1210019). TECZn.S / Valor: 1210019).

### For further information:

#### Tecan Group

Dr. Rudolf Eugster  
Chief Financial Officer

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

Martin Brändle  
Head of Corporate Communications &  
Investor Relations

Tel. +41 (0) 44 922 84 30  
Fax +41 (0) 44 922 88 89