

## **Tecan publishes 2009 Annual Report and notice of its Annual General Meeting**

**Männedorf, Switzerland, 30 March 2010** – Today the Tecan Group (SIX Swiss Exchange: TECN) announced the publication of its 2009 Annual Report on the company website [www.tecan.com](http://www.tecan.com) under Investor Relations. The Annual Report may also be requested in printed format. Tecan has already published its 2009 financial results in a press release and at the Press Briefing on Annual Results and Analysts' Conference held on 4 March 2010.

In addition, today the notice of the Tecan Group AG Annual General Meeting was published and dispatched. The Annual General Meeting will take place on 21 April 2010. This notice may also be found on the company website.

### **About Tecan**

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

### **For further information:**

#### **Tecan Group**

Martin Brändle

Head of Corporate Communications & Investor Relations

Tel. +41 (0) 44 922 84 30

Fax +41 (0) 44 922 88 89

[investor@tecan.com](mailto:investor@tecan.com)

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