# TECAN.

# **Press Release**

# Tecan publishes 2010 Annual Report and notice of its Annual General Meeting

**Männedorf, Switzerland, 29 March 2011** – Today the Tecan Group (SIX Swiss Exchange: TECN) announced the publication of its 2010 Annual Report on the company website <a href="www.tecan.com">www.tecan.com</a> under Investor Relations. The Annual Report may also be requested in printed format. Tecan has already published its 2010 financial results in a press release and at the Press Briefing on Annual Results and Analysts' Conference held on 10 March 2011.

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 19 April 2011. This notice may also be found on the company website.

#### **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 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 (OEM), 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 2010, Tecan generated sales of CHF 371 million (USD 356 million; EUR 269 million). Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TK: TECN/Reuters: TECZn.S/ ISIN: 12100191.

## 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
www.tecan.com